Agilent 1220 Infinity LC





© Agilent Technologies, Inc. 2010-2012

本マニュアルは米国著作権法およ び国際著作権法によって保護され ており、Agilent Technologies, Inc. の書面による事前の許可な く、本書の一部または全部を複製 することはいかなる形式や方法 (電子媒体による保存や読み出し、 外国語への翻訳なども含む)にお いても、禁止されています。

マニュアル番号

G4280-96016

エディション

05/2012

Printed in Germany

Agilent Technologies Hewlett-Packard-Strasse 8 76337 Waldbronn

本製品は、システムが適切な規制 機関で登録を受け関連する規制に 準拠している場合、ビトロ診断シ ステムのコンポーネントとして使 用できます。それ以外の場合は、 一般的な実験用途でのみ使用でき ます。

保証

このマニュアルに含まれる内容は 「現状のまま」提供されるもので、 将来のエディションにおいて予告 なく変更されることがあります。 また、Agilent は、適用される法 律によって最大限に許可される範 囲において、このマニュアルおよ びそれに含まれる情報に関して、 商品性および特定の目的に対する 適合性の暗黙の保証を含みそれに 限定されないすべての保証を明示 的か暗黙的かを問わず一切いたし ません。Agilent は、このマニュ アルまたはそれに含まれる情報の 所有、使用、または実行に付随す る過誤、または偶然的または間接 的な損害に対する責任を一切負わ ないものとします。Agilent とお 客様の間に書面による別の契約が あり、このマニュアルの内容に対 する保証条項がこの文書の条項と 矛盾する場合は、別の契約の保証 条項が適用されます。

技術ライセンス

このマニュアルで説明されている ハードウェアおよびソフトウェア はライセンスに基づいて提供さ れ、そのライセンスの条項に従っ て使用またはコピーできます。

安全に関する注意

注意

注意は、危険を表します。これは、正しく実行しなかったり、指示を順守しないと、製品の損害または重要なデータの損失にいたるおそれがある操作手順や行為に対する注条 やを十分に理解し、条件が視たされるまで、注意を無視して先に進んではなりません。

警告

警告は、危険を表します。こ れは、正しく実行しなかった り、指示を順守しないと、人 身への傷害または死亡にいた るおそれがある操作手順や行 為に対する注意を喚起します。 指示された条件を十分に理解 し、条件が満たされるまで、 警告を無視して先に進んでは なりません。



本書の内容

本書では、Agilent 1220 Infinity LC システムコンフィグレーションにつ いて説明します

- G4286B
- G4288B/C
- G4290B/C
- G4294B

1 概要

この章では、Agilent 1220 Infinity LC の使用可能なコンフィグレーションの概要を説明します。

2 設置要件と仕様

この章では、環境要件、物理的仕様、そして性能仕様について説明します。

3 据付

この章では、出荷内容と設置の概要を示します。

5 溶媒送液システムの説明

この章では、溶媒送液システム(ポンプとオプションのデガッサ)の操作 原理の概要を説明します。

6 注入システムの説明

この章では、注入システム(マニュアルインジェクタおよびオート サンプ ラ)の操作原理の概要を説明します。

7 カラムオーブンの説明

この章では、カラムオーブンの操作原理の概要を説明します。

本書の内容

8 検出器の説明

この章では、検出器の操作原理の概要を説明します。

9 テスト機能とキャリブレーション

この章では、Instrument Utilities ソフトウェアまたは Lab Advisor で 利用できるテスト、キャリブレーションおよびツールについて説明します。

10 エラー情報

この章では、表示されるエラーメッセージに関する情報を提供し、考えられる原因とそれらを解決するための提案を示します。

11 メンテナンス

この章では、機器のメンテナンスに関する一般情報を説明します。

12 メンテナンス用部品

この章では、メンテナンス用部品について説明します。

13 Agilent 1220 Infinity LC のアップグレード

この章では、LC システムのアップグレード用の情報を説明します。

14 ケーブルの識別

この章では、1200 シリーズの HPLC モジュールに使用されるケーブルにつ いて説明します。

1 概要 9

```
Agilent 1220 Infinity LC のコンフィグレーション 10
Agilent 1220 Infinity LC VL コンフィグレーション 11
EMF (Early Maintenance Feedback) 12
```

2 設置要件と仕様 15

設置について 16 物理的仕様 19 性能仕様 20

3 据付 31

システムの開梱 32 ハードウェアの据付方法 37 機器のクロマトデータシステムへの接続とコンフィグレーショ ン 46 Agilent 1220 Infinity LC と PC の接続 47 機器ユーティリティ / LabAdvisor ソフトウェア 49 アップグレード インストール後の機器の コンフィグレーショ ン 50 システムのプライミングと「据え付けチェックリスト」の実行 51 「チェックアウト分析」の実行 52

4 LAN コンフィグレーション 53

最初の必要事項 54 TCP/IP パラメータ コンフィグレーション 56 コンフィグレーション スイッチ 57 初期化モード選択 58 動的ホストコンフィグレーションプロトコル(DHCP) 62 リンクコンフィグレーション選択 66 BootP を使用した自動コンフィグレーション 67 Bootp を使用した設定の永久保存 78 マニュアルコンフィグレーション 79

目次

- 5 溶媒送液システムの説明 85
 概要 86
 デガッサ 87
 操作原理 88
 圧縮率補正 92
 可変ストローク容量 94
 ポンプの使用 95
- 6 注入システムの説明 97
 マニュアルインジェクタ 98
 オート サンプラ 102
- 7 カラムオーブンの説明 117 カラムオーブン 118

8 検出器の説明 119

検出器タイプ 120 Agilent 1220 Infinity LC 可変波長検出器 (VWD) 121 Agilent 1220 Infinity ダイオードアレイ検出器 (DAD) 122 適切なカラムとフローセルの組合せ 140

9 テスト機能とキャリブレーション 145
Agilent 1220 Infinity LC システム 147
溶媒送液システム 149
オート サンプラ 160
カラムオーブン 168
可変波長検出器 (VWD) 170
ダイオードアレイ検出器 (DAD) 180

- 10 エラー情報 205
 - エラー メッセージ内容 208 一般エラー メッセージ 209 ポンプのエラー メッセージ 219 オート サンプラのエラー メッセージ 237 一般検出器エラー メッセージ 252 VWD 検出器エラーメッセージ 257 DAD 検出器エラー メッセージ 267

11 メンテナンス 277

PM の作業範囲とチェックリスト 279 注意と警告 280 溶媒送液システム 282 マニュアルインジェクタ 303 オート サンプラ 307 可変波長検出器 (VWD) 327 ダイオードアレイ検出器(DAD) 338 HPLC システムでの藻の繁殖 359 モジュールのファームウェアの交換 361

12 メンテナンス用部品 363

1220 Infinity LC システム364溶媒送液システム366注入システム375カラムオーブン383検出器384

 13 Agilent 1220 Infinity LC のアップグレード
 391

 オーブンのアップグレード
 392

14 ケーブルの識別 393 ケーブル概要 394 アナログケーブル 396 リモートケーブル 398 BCD ケーブル 402 CAN/LAN ケーブル 405 Agilent 1200 モジュールから PC 406

15 付録 407

安全に関する一般的な情報 408 溶媒情報 411 無線干渉 413 紫外線放射 414 騒音レベル 415 廃電気電子機器(WEEE)指令(2002/96/EC) 416 ホルミウムフィルタの波長の証明書 417 Agilent Technologies のインターネットサービス 418



1220 Infinity LC

概要

Agilent 1220 Infinity LC のコンフィグレーション10Agilent 1220 Infinity LC VL コンフィグレーション11EMF (Early Maintenance Feedback)12ポンプの EMF カウンタ12オート サンプラの EMF カウンタ13可変波長型検出器の EMF カウンタ14ダイオードアレイ検出器の EMF カウンタ14

この章では、Agilent 1220 Infinity LC の使用可能なコンフィグ レーションの概要を説明します。



1 概要 Agilent 1220 Infinity LC のコンフィグレーション

Agilent 1220 Infinity LC のコンフィグレーション

Agilent 1220 Infinity LC の使用可能なコンフィグレーション

Agilent 1220 Infinity LC では、4 つの異なるコンフィグレーションで使 用できます。使用可能なコンポーネントは、アイソクラティック ポンプ、 デュアル チャンネル グラジエントポンプ (デガッサ付き)、マニュアルイ ンジェクタ、オート サンプラ、カラムオーブン、検出器などです。各コン フィグレーションには、少なくともポンプ 1 台、注入システム 1 台、検 出器 1 台が含まれ、Agilent Instrument Utilities ソフトウェアを搭載 しています。



溶媒切り替えバルブ(SSV)アップグレードキット(G4280-68708)が使用 可能です。

Agilent 1220 Infinity LC VL コンフィグレーション

Agilent 1220 Infinity LC VL の使用可能なコンフィグレーション

Agilent 1220 Infinity LC VL は、2 つの異なるコンフィグレーションで 使用できます。使用可能なコンポーネントは、アイソクラティック ポン プ、デュアル チャンネル グラジエントポンプ (デガッサ付き)、マニュア ルインジェクタ、オート サンプラ、カラムオーブン、検出器などです。各 コンフィグレーションには、少なくともポンプ 1 台、注入システム 1 台、 検出器 1 台が含まれ、Agilent Instrument Utilities ソフトウェアを搭 載しています。



溶媒切り替えバルブ(SSV)アップグレードキット(G4280-68708)が使用 可能です。

1

EMF (Early Maintenance Feedback)

ポンプの EMF カウンタ

EMF カウンタの EMF リミットはユーザーが設定可能なため、必要に応じて EMF (Early Maintenance Feedback) 機能を調整できます。ポンプコンポー ネントの摩耗は分析条件によって異なります。そのため、機器の固有運転 条件に基づき、上限値の定義を決定する必要があります。

Agilent 1220 Infinity LC ポンプは、ポンプヘッド用の一連の EMF カウ ンタを装備しています。各カウンタは、そのポンプが使用されるたびに増 加します。カウンタの上限値を割り当てておき、その制限を超えた時点で ユーザー インターフェースにフィードバックすることができます。メンテ ナンスの終了後、各カウンタをゼロにリセットできます。ポンプは、以下 の EMF カウンタを装備しています。

ポンプ送液量カウンタ

ポンプ送液量カウンタには、カウンタの最後のリセット以降ポンプヘッド が送液した溶媒の総量が表示されます。ポンプ送液量カウンタに、EMF (最大)限度を割り当てることができます。限界値を超えると、ユーザーイ ンターフェイスに EMF フラグが表示されます。

シール摩耗カウンタ

シール摩耗カウンタには、圧力と流量(共にシール摩耗の原因となる)か ら得られた値が表示されます。この値は、カウンタがシールメンテナンス 後にリセットされるまで、ポンプの利用と共に増加します。両方のシール 摩耗カウンタに、EMF(最大)限度を割り当てることができます。限界値 を超えると、ユーザーインターフェイスに EMF フラグが表示されます。

概要 EMF (Early Maintenance Feedback)

1

オート サンプラの EMF カウンタ

EMF カウンタの EMF リミットはユーザーが設定可能なため、必要に応じて EMF (Early Maintenance Feedback)機能を調整できます。オート サンプ ラ コンポーネントの摩耗は分析条件によって異なります。そのため、機器 の固有運転条件に基づき、上限値の定義を決定する必要があります。

オート サンプラは、2 つの EMF カウンタを装備しています。各カウンタ は、オート サンプラが使用されるたびに増加します。カウンタの上限値を 割り当ておき、その制限を超えた時点でユーザー インターフェースに フィードバックすることができます。メンテナンスの終了後、各カウンタ をゼロにリセットできます。オート サンプラは、次の EMF カウンタを装 備しています。

注入バルブカウンタ

このカウンタには、最後にカウンタがリセットされてからの、インジェク ションバルブの切り換え総数が表示されます。

ニードル移動カウンタ

このカウンタには、最後にカウンタがリセットされてからの、シートへの ニードルの動作総数が表示されます。

可変波長型検出器の EMF カウンタ

EMF カウンタの EMF リミットはユーザーが設定可能なため、必要性に合わ せて EMF (Early Maintenance Feedback)機能を調整できます。ランプの 有効点灯時間は、分析条件(高感度分析、低感度分析、波長など)によっ て異なります。そのため、機器の固有運転条件に基づき、上限値の定義を 決定する必要があります。

検出器モジュールは、ランプ用の EMF カウンタを装備しています。カウン タは、ランプが使用されるたびに増加します。カウンタの上限値を割り当 てておき、その制限を超えた時点でユーザインタフェースにフィードバッ クすることができます。ランプの交換後、カウンタをゼロにリセットする ことができます。本検出器は、以下の EMF カウンタを装備しています。

重水素ランプ点灯時間

このカウンタには、重水素ランプの合計点灯時間が時間単位で表示されま す。

ダイオードアレイ検出器の EMF カウンタ

EMF カウンタの使用

EMF カウンタの EMF リミットはユーザーが設定可能なため、必要に応じて EMF 機能を調整できます。有効なメンテナンスサイクルは使用要件によっ て異なります。そのため、機器に固有の動作条件に基づいて最大リミット 値の定義を決定する必要があります。

EMF リミットの設定

EMF リミットの設定は、1 回または 2 回以上のメンテナンスサイクルにわ たって最適化します。最初にデフォルトの EMF リミット値を設定する必要 があります。性能の低下によってメンテナンスが必要であることがわかっ た場合は、EMF カウンタの表示値を書き留めておいてください。これらの 値(または表示された値より多少小さい値)を EMF リミットとして入力 し、EMF カウンタをゼロにリセットします。次に EMF カウンタがこの EMF リミットを超えると、EMF フラグが表示され、メンテナンスが必要な時期 であることを知らせます。



1220 Infinity LC

2 設置要件と仕様

設置について	•	16
電源につい	いて	16
電源コー	ド	17
設置スペー	ース	17
環境	18	
物理的仕様	19	
性能仕様	20	
仕様条件	29	9

この章では、環境要件、物理的仕様、そして性能仕様について説明 します。





設置について

機器を最適な性能で動作させるためには、適切な環境に設置することが重 要です。

電源について

Agilent 1220 Infinity LC のパワー サプライは、広範囲の入力電力に対応しています。したがって、機器には電圧スイッチはありません。

警告オフの状態でも、機器は部分的に通電しています。

フロント パネルの電源スイッチがオフになっている場合でも、パ ワー サプライは少量の電力を使用します。検出器の修理作業により 人身障害に至る恐れがあります。たとえば、検出器カバーが開いて いて機器が電源に接続されている場合の感電の危険などです。

→検出器を電源供給ラインから切り離すには、電源コードを抜いて ください。

- 警告
 機器への不適正な入力電圧
 仕様より高い入力電圧に接続すると、感電の危険が生じたり、機器
 を損傷することがあります。
 → 使用する機器は、指定された入力電圧だけに接続してください。
 - 注意 緊急時に備えて、いつでも電源供給ラインから装置を切り離すことができ るようにしておく必要があります。

電源ケーブルは手の届きやすい場所に配置し、装置を入力電圧からすぐ切り離せるようにしてください。

→ ケーブルを抜けるように、機器の電源ソケットの後には十分な空間を確 保してください。

電源コード

機器には、オプションとして各種の電源コードが用意されています。どの 電源コードも、メス型側の形は同じです。電源コードのメス型を、機器の 左背面にある電源ケーブル コネクタに差し込みます。電源コードのオス型 側はコードによって異なり、各使用国または各地域のアース付きコンセン トに合わせて設計されています。

<u>警告</u> 接地不備または指定外の電源コードの使用

接地しなかったり、指定外の電源コードを使用すると、感電や回路 がショートすることがあります。

- → 接地していない電源を使用して機器を操作しないでください。
- → また、使用する地域に合わせて設計された Agilent Technologies の電源コード以外は使用しないでください。

■ Agilent が提供したものでないケーブルの使用

Agilent Technologies が提供したもの以外のケーブルを使用する と、電子コンポーネントが損傷したり、人体に危害が及ぶことがあ ります。

→ 安全基準または EMC 規格のコンプライアンスと正しい動作を確実 にするために、Agilent Technologies 製以外のケーブルは使用し ないでください。

設置スペース

Agilent 1220 Infinity LC は、機器をほとんどの机やラボの作業台に置く ことができるような寸法と重量になっています。通気と配線用に、本機器 の両側に 2.5 cm (1.0 in) および、背面に約 8 cm (3.1 in) の追加ス ペースが必要です。

Agilent 1220 Infinity LC を支えるための作業台が、機器の重量に耐える よう設計されているか確認します。

Agilent 1220 Infinity LC は直立状態で操作する必要があります。

警告

環境

注意

Agilent 1220 Infinity LC は、以下のセクションに記載されている周囲温 度および相対湿度の仕様範囲内で動作します。

ASTM ドリフトテストには、1 時間にわたる測定で温度変化が 2 °C/hour (3.6 °F/hour)未満になる環境条件が必要です。弊社が作成したドリフト 仕様は、上記の条件に基づいています。周囲温度変化が大きくなると、ド リフトも大きくなります。

ドリフトパフォーマンスは、温度変化をコントロールすることで改善できます。最高のパフォーマンスを実現するには、温度変化の頻度と幅を最小限に抑え、1°C/hour(1.8°F/hour)未満に保ちます。ただし、1分以内程度の短時間の変動は無視できます。

モジュール内の結露

結露によってシステムの電気回路が損傷することがあります。

- → 温度変化によってモジュール内に結露が発生する可能性がある環境条件 では、モジュールの保管、輸送、または使用を行わないでください。
- → 寒冷な天候下でモジュールが出荷された場合は、結露が発生しないよう に、オートサンプラを梱包箱に入れたままゆっくり室温まで温度を上げ てください。

物理的仕様

タイプ	仕様	コメント
重量	30 kg (66 lbs) G4294B: 43 kg (94 lbs)	
寸法 (高さ × 幅 × 奥行き)	$\begin{array}{l} 640 \ \times \ 370 \ \times \ 420 \ \mathrm{mm} \\ (25.2 \ \times \ 14.6 \ \times \ 16.5 \ \text{l}) \\ 64294 \\ \mathrm{B}: \\ 640 \ \times \ 370 \ \times \ 485 \ \mathrm{mm} \\ (25.2 \ \times \ 14.6 \ \times \ 19.1 \ \text{l}) \\ \end{array}$	
入力電圧	100 - 240 VAC, \pm 10 %	広範囲の電圧に対応
電源周波数	50 または 60 Hz、± 5 %	
消費電力	240 VA / 210 W / 717 BTU	最大値
使用周囲温度	4-55 ° C (39-131 ° F)	
保管周囲温度	-40 - 70 $^{\circ}$ C (-4 - 158 $^{\circ}$ F)	
湿度	$<$ 95 % r.h. \checkmark 40 $^{\circ}$ C (104 $^{\circ}$ F)	結露なし
使用高度	最高 2000 m (6562 ft)	
保管高度	最高 4600 m (15091 ft)	モジュールを保管で きる高度
安全規格: IEC、CSA、UL	設置クラス Ⅱ、汚染度 2	室内使用専用。

表 1 物理的仕様

性能仕様

Agilent 1220 Infinity LC の性能仕様

タイプ	仕様
安全機能	拡張診断機能、エラー検出と表示、 リーク検出、安全なリーク処理、ポ ンプ システムのシャットダウン用 リーク出力シグナル。主要なメンテ ナンス領域における低電圧。
コントロールおよびデータ評価	Agilent EZChrom Compact, Agilent ChemStation, Agilent Instrument Utilities, Agilent Lab Advisor
通信	コントローラエリアネットワーク (CAN)、RS232C、APG リモート:レ ディ、開始、終了、シャットダウン シグナル、LAN
GLP 機能	EMF (Early Maintenance Feedback)、 メンテナンスおよびエラーの電子記 録

表 2 Agilent 1220 Infinity LC の性能仕様

Agilent 1220 Infinity LC ポンプの性能仕様

表 3 Agilent 1220 Infinity LC ポンプの性能仕様

タイプ	仕様
ハイドロリック システム	弊社独自のサーボ制御方式可変ストロークドライブ、 フローティングプランジャ、パッシブインレットバ ルブを搭載したデュアルプランジャ直列型ポンプ
設定可能な流量範囲	0.001 - 10 mL/min (0.001 mL/min 単位)
流量範囲	0.2 - 10.0 mL/min
流量精度	一定室温でのリテンションタイムに基づき、≤0.07 % RSD、または< 0.02 min SD のいずれか大きい方
流量真度	± 1 % または 10 μL/min のいずれか大きい方、脱 気 H ₂ 0、80 - 100 bar、1 mL/min、一定室温
圧力	動作範囲 0 - 60 MPa (0 - 600 bar、0 - 8820 psi) 最高 5 mL/min 動作範囲 0 - 20 MPa (0 - 200 bar、0 - 2950 psi) 最高 10 mL/min
圧力の脈動	< 2 % 振幅(通常< 1.3 %)、1 mL/min イソプロパ ノール、すべての圧力> 1 MPa(10 bar)
圧縮率補正	移動相の圧縮率に応じて、切り換え可
推奨 pH 範囲	1.0 - 12.5 、pH < 2.3 の溶媒は、ステンレス ス チールを侵す酸を含まないこと
グラジエント形成 (オプション)	弊社独自の高速比例バルブを使用した低圧バイナリ 溶媒混合 / グラジエント機能
ディレイ ボリューム	600 - 900 μL、背圧によって異なる。測定時水量 1 mL/min (水/カフェイントレーサー)
組成範囲	0 - 95 % または 5 - 100 、切り換え可
組成精度	一定室温でのリテンションタイムに基づき、 < 0.2 % RSD または< 0.4 min SD のいずれか大き い方、1 mL/min

Agilent 1220 Infinity LC ポンプ VL の性能仕様

タイプ	仕様
ハイドロリック システム	弊社独自のサーボ制御方式可変ストロークドライブ、 フローティングプランジャ、パッシブインレットバ ルブを搭載したデュアルプランジャ直列型ポンプ
設定可能な流量範囲	0.001 - 10 mL/min、0.001 mL/min 単位
流量範囲	0.2 - 10 mL/min
流量精度	一定室温でのリテンションタイムに基づき、< 0.07 % RSD、または< 0.02 min SD のいずれか大き い方
流量真度	± 1 % または 10 μL/min のいずれか大きい方
圧力	動作範囲 0 - 40 MPa (0 - 400 bar、0 - 5880 psi) 最高 5 mL/min 動作範囲 0 - 20 MPa (0 - 200 bar、0 - 2950 psi) 最高 10 mL/min
圧力の脈動	< 2 % 振幅 (通常 < 1 %)、1 mL/min イソプロパ ノール、すべての圧力 > 1 MPa (10 bar)
圧縮率補正	移動相の圧縮率に応じて、切り換え可
推奨 pH 範囲	1.0 - 12.5 、pH < 2.3 の溶媒は、ステンレス ス チールを侵す酸を含まないこと
グラジエント形成 (オプション)	弊社独自の高速比例バルブを使用した低圧クォータ ナリ混合 / グラジエント機能、ディレイボリューム 800 ~ 1100 µL、背圧により異なる
組成範囲	0 ~ 95 % または 5 ~ 100 %、切り換え可
組成精度	< 0.2 % RSD、0.2 および 1 mL/min

表 4 Agilent 1220 Infinity LC ポンプ VL の性能仕様

Agilent 1220 Infinity LC オート サンプラの性能仕様

タイプ	仕様
圧力	動作範囲 0 - 60 MPa (0 - 600 bar、0 - 8820 psi)
注入範囲	0.1 - 100 μL0.1 μL 単位、マルチ注入で最高 1500 μL(ハードウェアの調整が必要)
繰り返し注入回数	バイアル 1 本で 1 ~ 99 回
精度	< 0.25 % RSD は下限 5 - 100 μL、< 1 % RSD 1 - 5 μL 容量可変
最小サンプル容量	1 μL 下限 5 μL サンプルで 100 μL マイクロバイ アル、または 1 μL、下限 10 μL サンプルで 300 μL マイクロバイアル
キャリーオーバ	通常< 0.1 %、外部ニードルのクリーニングで < 0.05 %
サンプルの粘性の範囲	0.2 - 50 cp
サンプルのキャパシ ティファクタ	1 トレイに 100 × 2 mL バイアル ½ トレイに 40 × 2 mL バイアル ½ トレイに 15 × 6 mL バイアル (Agilent バイ アルのみ)
注入サイクルタイム	通常 50 s 吸引速度と注入量により異なる

表 5 Agilent 1220 Infinity LC オート サンプラの性能仕様

Agilent 1220 Infinity LC カラムオーブンの性能仕様

タイプ	仕様
温度範囲	5 ° C 室温以上~ 60 ° C 5 ° C 室温以上~ 80 ° C (極小 FW Rev. B.06.50)
温度の安定性	± 0.15 ° C、組成および流速一定
温度の正確さ	\pm 0.8 $^{\circ}$ C
カラム容量	25 cm カラム 1
内部容量	6 µ L

表 6 Agilent 1220 Infinity LC カラムオーブンの性能仕様

Agilent 1220 Infinity LC VWD の性能仕様

表 7 Agilent 1220 Infinity LC VWD の性能仕様

タイプ	仕様	コメント
検出器タイプ	ダブルビーム分光光度計	
光源	重水素ランプ	
波長範囲	190 - 600 nm	
ノイズ	0.35・10 ⁻⁵ AU で 230 nm	時定数2秒、指定条件下
ドリフト	$3\cdot10^{-4}$ AU/hr で 254 nm	表下の注記を参照してくだ さい。
直線性	> 2 AU(5 %)上限値	表下の注記を参照してくだ さい。
波長真度	\pm 1 nm	重水素ラインによるセルフ キャリブレーションおよび 酸化ホルミウム フィルタに よるベリフィケーション。
最大データ速度	80 Hz	
バンド幅	6.5 nm 通常	
フローセル	標準: 14 µL 容量、10 mm セル光路長および 40 bar (588 psi) 最大圧力 高圧: 14 µL 容量、10 mm セル光路長および 400 bar (5880 psi) 最大圧力 セミマイクロ: 5 µL 容 量、6 mm セル光路長およ び 40 bar (588 psi) 最大圧力 マイクロ: 2 µL 容量、 3 mm セル光路長および 40 bar (588 psi) 最大圧力	コンポーネントレベルの修 理が可能です。



注記

ASTM: 『液体クロマトグラフィに使用する可変波長型光度検出器の実施基準』 基準条件:セル光路長 10 mm、レスポンス タイム 2 s、流量 1 mL/min LC ク ラスのメタノール。直線性は、カフェインで 272 nm にて測定。

Agilent 1220 Infinity LC DAD の性能仕様

タイプ	仕様	コメント
検出器タイプ	1024 素子ダイオードアレイ	
光源	重水素ランプとタングステ ンランプ	UV ランプには、ランプの 標準的な情報を含む RFID (無線認識)タグが取り 付けられています。
シグナル数	8	
最大サンプリング スピード	80 Hz	
短期ノイズ (ASTM) シングル およびマルチ 波長	< ± 0.7・10 ⁻⁵ AU、 254/4 nm および 750 nm、 TC 2 s	後の「仕様条件」を参照 してください。
ドリフト	$<~0.9\cdot~10^{\rm -3}$ AU/h, 254 nm	後の「仕様条件」を参照 してください。
線形吸光度範囲	> 2 AU (5 %), 265 nm	後の「仕様条件」を参照 してください。
波長範囲	190 - 950 nm	
波長真度	\pm 1 nm	重水素ラインによるセル フキャリブレーションお よび酸化ホルミウム フィ ルタによるベリフィケー ション。
スリット幅	1、2、4、8、16 nm	プログラム式スリット
ダイオード幅	< 1 nm	

表 8 性能仕様

表	8	性能	仕様
~	-	1-1-1-1-1-1	

タイプ	仕様	コメント
フローセル	標準: 13 µL 容量、10 mm セル光路長および 120 bar (1740 psi) 圧力最大	フローセルには、セルの 標準的な情報を含む RFID (無線認識) タグが取り 付けられています。 pH 範囲 1.0 ~ 9.5
時間プログラム可 能	波長、極性、ピーク幅、ラ ンプ帯域幅、自動バランス、 波長範囲、スレッショルド、 スペクトル保存モード	

仕様条件

ASTM: 『液体クロマトグラフィに使用する可変波長型光度検出器の実施基準』

基準条件:セル光路長 10 mm、波長 254 および 750 nm、リファレンス波 長 360 nm/100 nm 使用、スリット幅 4 nm、時定数 2 s (レスポンス タイ ム 4 s に相当)、流量 1 mL/min LC クラスのメタノール。

直線性: 直線性は、カフェインで 265 nm/4 nm、スリット幅 4 nm および TC 2 s (または RT 4 s)、光路長 10 mm を使用して測定。

環境条件については、「環境」を参照してください。

注記 この仕様は標準 RFID タグ付きランプ(2140-0820)を基準にしているので、 その他のランプタイプを使用する場合は実現できない可能性があります。

注記 機器の近くでモバイルデバイスを使用すると、検出器の短期ノイズレベルに影響することがあります。

ASTM ドリフトテストには、1 時間にわたる測定で温度変化が 2 ° C/hour (3.6 ° F/hour)未満であることが必要です。弊社が作成したドリフト仕様 は、上記の条件に基づいています。周囲温度変化が大きくなると、ドリフ トも大きくなります。ドリフトパフォーマンスは、温度変化をコントロー ルすることで改善できます。最高のパフォーマンスを実現するには、温度 変化の頻度と幅を最小限に抑え、1 ° C/hour (1.8 ° F/hour)未満に保ち ます。ただし、1 分以内程度の短時間の変動は無視できます。

パフォーマンステストは、完全にウォームアップした光学ユニット(> 2 時間)で行う必要があります。ASTM 測定では、テスト開始の少なくとも 24 h 前に検出器の電源を入れる必要があります。

時定数 / レスポンスタイム

ASTM E1657-98 『液体クロマトグラフィに使用するの可変波長型光度検出 器テストの実施基準』に従い、係数 2.2 を掛けることで、時定数をレスポ ンスタイムに変換します。





1220 Infinity LC

-据付

3

システムの開梱 32 梱包明細リスト 32 ハードウェアの据付方法 37 据付方法 37 Agilent 1220 Infinity LC の設置 39 1220 Infinity LC の接続の識別 44 機器のクロマトデータシステムへの接続とコンフィグレーショ ン 46 Agilent 1220 Infinity LC と PC の接続 47 機器ユーティリティ / LabAdvisor ソフトウェア 49 アップグレード インストール後の機器の コンフィグレーショ ン 50 システムのプライミングと「据え付けチェックリスト」の実 行 51 「チェックアウト分析」の実行 52

この章では、出荷内容と設置の概要を示します。

注記

Agilent 1220 Infinity LC システムを設置するため、インストールガイドに 段階的に従うことを強くお勧めします。



3 据付 システムの開梱

システムの開梱

梱包材の外部に損傷の形跡がある場合は、Agilent Technologies の営業所 およびサービスオフィスまで即ご連絡ください。サービス担当者に、 Agilent 1220 Infinity LC が配送中に損傷を受けた可能性があることをご 通知ください。

注意 損傷の形跡

→ Agilent 1220 Infinity LC を取り付けようとしないでください。

梱包明細リスト

梱包明細リスト

Agilent 1220 Infinity LC と一緒にすべての部品と器材が納品されたこと を確認してください。梱包明細リストを以下に示します。不足品または破 損品があった場合は、Agilent Technologies の営業所およびサービスオ フィスまでご連絡ください。

表 9 Agilent 1220 Infinity チェックリスト

説明	個数
Agilent 1220 Infinity LC	1
電源ケーブル	1
フローセル	据付済み
機器ユーティリティ DVD	1
インストールガイド	1
アクセサリキット (下記参照)	1

G4286B 用アクセサリキット内容

部品番号	説明
G4286-68755	アクセサリキット一式
0100-2562	フィッティング、一体型、指締め式
0890-1195	PTFE チューブ、内径 0.052 in
0890-1711	可撓管 (排出用)、3 m
5023-0203	クロスオーバーネットワークケーブル、シールド付き、 3 m (ポイントツーポイント接続用)
5062-8535	廃液アクセサリキット
5188-2758	PTFE/シリコンセプタム、16 mm、プレスリット、 100 個入り (納品個数 0.010)
5190-1501	シリンジ、50.0 μL、FN、LC チップ
9301-0411	シリンジ、プラスチック製
9301-1337	シリンジアダプタ
9301-1377	スクリューキャップバイアル、透明、6 mL 100 個入り (納品個数 0.010)
9301-1379	スクリューキャップ、6 mL バイアル用、100 個入り (納品個数 0.010)
9301-1420	溶媒ボトル、透明
G1311-60003	ボトルヘッドアセンブリ

G4288B/C 用アクセサリキット内容

部品番号	説明
G4288-68755	アクセサリキット一式
0100-2562	フィッティング、一体型、指締め式
0890-1195	PTFE チューブ、内径 0.052 in
0890-1711	可撓管 (排出用)、3 m
5023-0203	クロスオーバーネットワークケーブル、シールド付 き、3 m(ポイントツーポイント接続用)
5062-8535	廃液アクセサリキット
5188-2758	PTFE/シリコンセプタム、16 mm、プレスリット、100 個入り (納品個数 0.010)
5190-1501	シリンジ、50.0 µL、FN、LC チップ
9301-0411	シリンジ、プラスチック製
9301-1337	シリンジアダプタ
9301-1377	スクリューキャップバイアル、透明、6 mL 100 個入り (納品個数 0.010)
9301-1379	スクリューキャップ、6 mL バイアル用、100 個入り (納品個数 0.010)
9301-1420	溶媒ボトル、透明
9301-1450	溶媒ボトル、褐色
G1311-60003 (2x)	ボトルヘッドアセンブリ

G4290B/C、G4294B 用アクセサリキット内容

部品番号	説明
G4290-68755	アクセサリキット一式
0100-2562	フィッティング、一体型、指締め式
0890-1195	PTFE チューブ、内径 0.052 in
0890-1711	可撓管 (排出用)、3 m
5023-0203	クロスオーバーネットワークケーブル、シールド付 き、3 m(ポイントツーポイント接続用)
5062-8535	廃液アクセサリキット
9301-0411	シリンジ、プラスチック製
9301-1337	シリンジアダプタ
9301-1420	溶媒ボトル、透明
9301-1450	溶媒ボトル、褐色
G1311-60003 (2x)	ボトルヘッドアセンブリ

Agilent 1220 Infinity LC 用ツールキットオプション

部品番号	説明
G4296-68715	ツールキットー式
0100-1710	チューブ接続用取り付けツール
8710-0510 (2x)	1/4 インチ × 5/16 インチのレンチ
8710-1924	両ロスパナ 14 mm
8720-0025	スパナ、1/2 inch & 9/16 inch
01018-23702	インサートツール
8710-2392	六角レンチ4 mm15 cm 、T 字型ハンドル
8710-2394	六角レンチ、9/64 インチ15 cm 長、T 字型ハンドル
8710-2411	六角レンチ3 mm12 cm 長
8710-2412	六角レンチ2.5 mm、15 cm 長、ストレートハンドル付き
8710-0899	Pozidriv ドライバ
ハードウェアの据付方法

ハードウェアの据付方法

据付方法

標準据付方法

(据付中は、モジュールにオプションのハード ウェアは追加されません。) アップグレードキットの設置を含めた据付方法

(オーブンのアップグレードキット/マニュアル インジェクタから ALS へのアップグレードキッ ト/アイソクラティックからグラジエントへの アップグレードキット)

注記

グラジエントシステムのアップグレードキット および ALS アップグレードキットのインストー ルは、Agilent の訓練を受けたサービス担当の みが行います。

ハードウェアの据付方法

標準据付方法



アップグレードキットの設置を含めた据付方法



注記

溶媒切り替えバルブ (SSV) オプションの据付に は、新しい機器タイプのコンフィグレーション は必要ありません。CDS のみ、SSV の設定が必 要です。

ハードウェアの据付方法

Agilent 1220 Infinity LC の設置

- 1 開梱し、内容を明細リストと比較してすべて揃っていることを確認 します。
- 2 Agilent 1220 Infinity LC を作業台の上に設置します。
- 3 リリース ボタン (両側) を押して、両方のフロント カバー (上下) を 取り外します。



図 1 フロント カバーのメカニズム

ハードウェアの据付方法



4 2 個の配送用発泡材を取り外します。

図 2 配送用発泡材の取り外し

『41 ページ 図 3』は、フロント カバーを取り外した、フル装備の 1220 Infinity LC システムの内容を示します。 (表示されているモジュール のタイプは、G4290B)

居付 3

ハードウェアの据付方法



図 3 1220 Infinity LC システムの概要

注記

追加のオプションまたはアップグレードキットは、すべての溶媒の流路を設置 する前に取り付けます。Agilent 1220 Infinity LC モジュール情報のコンフィ グレーションについては、『「アップグレード インストール後の機器の コン フィグレーション」50ページ』に記載されています。

オプションとアップグレードキットの据付方法の詳細については、Agilent 1220 Infinity LC モジュールのマニュアルを参照してください。

5 0.5 L の HPLC クラスの水を入れた溶媒ボトルを溶媒トレイに設置 します。

ハードウェアの据付方法

6 溶媒インレットフィルタを溶媒ボトルのボトル ヘッド アセンブリの端 に設置します(下図を参照)。



図 4 ボトル ヘッド アセンブリおよび溶媒ボトル

- 7 チューブが水でいっぱいになるまで、シリンジ (9301-044) および シリ ンジアダプタ (9301-1337) (アクセサリキットの一部)を使用して水を 入れます。
- 8 ボトル ヘッド アセンブリのコネクタを以下に接続します(『42ページ図4』項目1と2を参照)。
 - パッシブインレットバルブ (アイソクラティックポンプ)
 - デガッサ注入口チャンネル A (グラジエントポンプ)
- 9 廃液チューブとフローセルアウトレットに付属するフィッティング (アクセサリキットの一部)を接続し、もう一方の端を適切な溶媒廃液 コンテナに接続します(『42ページ図5』を参照)。



図 5 VWD モジュールの接続

ハードウェアの据付方法

- 10 波形廃液チューブ(アクセサリキットの一部)を VWD リークトレイ出口 のアダプタに取り付け、適切な廃液コンテナに導きます(『42 ページ 図 5』を参照)。
- 11 廃液チューブ(アクセサリキットの一部)をパージ バルブ出口のアダ プタに接続し、もう一方の端を廃液コンテナに接続します。
- **12** Agilent 1220 Infinity LC と PC 間のネットワークを接続します。

注記 機器のネットワーク接続を確立する方法の詳細については、以下を参照してく ださい。『「Agilent 1220 Infinity LC と PC の接続」47 ページ』または 『「LAN コンフィグレーション」53 ページ』

- **13** モジュールの前部にある電源プッシュボタンが離れていることを確認 します(『42ページ図5』を参照)。ここで、電源コードを Agilent 1220 Infinity LC と電源供給ラインに接続します。
- 14 モジュールの電源を入れる前に、発泡材をすべて取り外したことを確認 します(『40 ページ図2』を参照)。次に電源用プッシュボタンでモ ジュールの電源を入れます。

1220 Infinity LC の接続の識別

Agilent 1220 Infinity LC と VWD

『44 ページ 図 6』は、1220 Infinity LC 機器と VWD 間の可能な接続の概 要を示します。



図 6 Agilent 1220 Infinity LC と VWD の接続

ハードウェアの据付方法

Agilent 1220 Infinity LC と DAD

『45 ページ 図 7』は、1220 Infinity LC 機器と DAD の可能な接続の概要 を示します。G4294B では、DAD メインボード上のコンフィグレーション ディップ スイッチを使用する必要があります。この場合、それが通信ホス トとなるからです。短い CAN ケーブルは、DAD と他の 1220 モジュール間 の通信接続用です。



図 7 Agilent 1220 Infinity LC と DAD の接続

機器のクロマトデータシステムへの接続とコンフィグレーション

機器のクロマトデータシステムへの接続とコンフィグ レーション

- 1 クロマトデータシステム (CDS) をインストールします。CDS に付属のイ ンストールマニュアルを参照してください。
- **2** CDS を開始します。
- **3** 機器コンフィグレーションの画面で、機器名(自由選択)と機器タイプ を入力します (Agilent Compact LC)。
- 4 モジュールのコンフィグレーションには、自動コンフィグレーションを 選択します。

Agilent 1220 Infinity LC と PC の接続

AGILENT 1220 Infinity LC は、デフォルトのネットワークコンフィグレー ション設定で工場から出荷されます。(コンフィグレーション ディップ スイッチ 7 & 8 は ON に設定されています)。これにより、クロスオー バーパッチケーブル(アクセサリキットの一部)を通じて、PC に速やかに 接続できます。

工場出荷時のデフォルトの IP アドレス: 192.168.254.11



注記

G4294B では、DAD 拡張ボードのコンフィグレーション ディップ スイッチを 使用して、LAN 接続を設定します。

Agilent 1220 Infinity LC と PC の接続

このデフォルト アドレスを使用して機器を PC に接続するためには、以下のように PC ネットワークを設定します。

IP:	192. 168. 254. 10
サブネットマスク:	255. 255. 255. 0
デフォルトゲートウェイ:	該当なし

2 Agilent 1220 Infinity LC の LAN-ポート (『47 ページ 図 8』) と PC の ネットワーク コネクタを、クロスオーバーパッチケーブルで接続しま す。

注記 クロスオーバーケーブルは、モジュールと PC 間の直接接続専用です。 Agilent 1220 Infinity LC をハブを通じてネットワークに接続する場合は、 最寄りのネットワーク管理者に相談してください。

機器をネットワークに接続する場合は、有効なネットワークアドレスを提 供できる最寄りのネットワーク管理者に相談することを強くお勧めします。 Agilent 1220 Infinity LC の LAN コンフィグレーションの詳細について は、1220 Infinity LC のユーザーマニュアルの「Installation」の章にあ る「LAN Configuration」を参照してください。

以下の詳細が記載されています。

- TCP/IP パラメータ コンフィグレーション
- ・ コンフィグレーション スイッチ
- 固定 IP アドレスの使用
- 個別の IP アドレスの設定方法

機器ユーティリティ / LabAdvisor ソフトウェア

機器ユーティリティ / LabAdvisor ソフトウェア

Agilent 1220 Infinity LC インストールプロセス中は、このソフトウェア を使用して、システムのフラッシングとシステムインストールチェックを 実行します(『「システムのプライミングと「据え付けチェックリスト」の 実行」51 ページ』を参照)。

- 1 ソフトウェア CD のインストールの手順に従い、機器ユーティリティま たは Lab Advisor ソフトウェアをインストールします。
- 2 機器をソフトウェアに設定し、それを接続します。

注記 ハードウェアのアップグレード (アイソクラティックからグラジエント、カラ ムオーブン、オート サンプラへのアップグレード)をするときは常に、機器 を機器ユーティリティ / Lab Advisor ソフトウェアで再設定する必要がありま す。そのため、続行する前に、『「アップグレード インストール後の機器の コ ンフィグレーション」50ページ』に記載されている手順に従ってください。

アップグレード インストール後の機器の コンフィグレーション

アップグレード インストール後の機器の コンフィグレーション

このステップは、以下のハードウェアのアップグレードキットを機器にインストールした場合のみ必要です。

- G4297A 1220 Infinity オーブンキット
- G4298A 1220 Infinity アップグレード マニュアルインジェクタから ALS
- G4299A 1220 Infinity アップグレード アイソクラティックからグラ ジェントポンプ
- 1 機器を機器ユーティリティ / Lab Advisor ソフトウェアへ接続します。
- 2 適用したハードウェアの変更内容に従って機器を設定します。 ソフトウェアリビジョン B.01.04 以下:
 - ツール > モジュール サービスセンター (任意のモジュール)

ソフトウェアリビジョン B.02.01 以上:

機器コントロール > コントロール(任意のモジュール) > コンバー
 ト デバイス タイプ

(たとえば、G4297A - 1220 Infinity オーブンキットを設置した場合は、 [オーブンを追加]を押します。

- **3** ソフトウェアから切り離し、機器を再起動します。
- 4 機器を機器ユーティリティ / Lab Advisor ソフトウェアへ再接続します。
- 5 ここで、据え付けチェックリスト(『「システムのプライミングと「据え 付けチェックリスト」の実行」51ページ』に記載)を実行します。
- 6 クロマトグラフ データシステムを開始するときは、自動コンフィグ レーションを使用して、新しいハードウェアのコンフィグレーションを CDS に設定します。

システムのプライミングと「据え付けチェックリスト」の実行

システムのプライミングと「据え付けチェックリスト」の 実行

以下に記載されるステップは、機器ユーティリティ / Lab Advisor ソフト ウェアを使用して実行します。

- すべてのチャンネルを HPLC クラスの水に接続し、パージポンプを使用 して溶媒チャンネルをフラッシングします。 ソフトウェアリビジョン B.01.04 以下:
 - ツール > ポンプ > パージポンプ
 - ソフトウェアリビジョン B.02.01 以上:
 - サービス & 診断 > ポンプ(ツールのチェックは必須) > パージポン
 プ
- すべての接続したチャンネルを、バブルフリーになるまで十分プライミングします。
- **3 機器コントロール**の機能性を使用して、システムを HPLC クラスの水で プライミングし、システムから空気を除去します。
 - 以下の条件を適用します。
 - パージ バルブ:閉じる
 - 流量: 2 mL/min
 - 時間: 5 min/channel
 - 設定ストローク: 100 µL
- 4 [サービスと診断] メニューの [据え付けチェックリスト] を実行し、 テスト結果を印刷します。
- 5 ステータスレポートを作成し印刷します。

「チェックアウト分析」の実行

「チェックアウト分析」の実行

- 1 クロマトデータシステムを開始します。
- 2 以下のパラメータでチェックアウトメソッドを作成します。
 - 流量:1 mL/min
 - 注入 容量: 20 µL (オート サンプラ)
 - オーブン温度:コントロールなし
 - VWD の波長: 254 nm
 - 分析時間: 1 min

このチェックアウト分析は、工場で据付済みの「リストリクション キャピラリ」を使用して実行します。

- 3 1ml の標準試料(たとえば、アセトン)を準備し、オート サンプラ ト レイのバイアルポジション 1 に設置します。 マニュアルインジェクタ コンフィグレーションには、20 µL の標準試料 をループに充填します。注入ループを少なくとも 3 倍満杯にします。
 - (例、少なくとも 60 µL を 20 µL サンプル ループで注入)
- 4 シングル ランを開始します。結果、シングルピークが確認されるはずです。
- 5 レポートを印刷します。
- 6 作成印刷したすべてのレポートをバインダーに保存します。 これで、Agilent 1220 Infinity LC のインストールが完了しました。



1220 Infinity LC

4

・ LAN コンフィグレーション

最初の必要事項 54 TCP/IP パラメータ コンフィグレーション 56 コンフィグレーション スイッチ 57 初期化モード選択 58 動的ホストコンフィグレーションプロトコル (DHCP) 62 一般情報 (DHCP) 62 セットアップ (DHCP) 64 リンクコンフィグレーション選択 66 BootP を使用した自動コンフィグレーション 67 Agilent BootP サービスについて 67 BootP サービスの使用 68 状況: LAN 通信を確立できない場合 68 BootP サービスのインストール 69 MAC アドレスを決定する 2 つの方法 71 Agilent BootP サービスによる IP アドレスの割り当 T 73 Agilent BootP サービスによる機器の IP アドレスの変 更 76 Bootp を使用した設定の永久保存 78 マニュアルコンフィグレーション 79 Telnet の使用 80



4 LAN コンフィグレーション 最初の必要事項

最初の必要事項

Agilent 1220 Infinity LC には、オンボードの LAN 通信インタフェース が装備されています。

今後の参照用に、MAC (メディアアクセス コントロール) アドレスを書き留めてください。LAN インタフェースの MAC アドレスまたはハードウェアアドレスは、世界中で唯一の ID です。別のネットワーク機器が同じハードウェアアドレスを持つことはありません。コンフィグレーション スイッチの隣の機器背面左側に MAC アドレスのラベルがあります。



検出器メインボードの部品番号 リビジョンコード、製造供給元、 アセンブリの年と週 MAC アドレス 生産国

図 9 MAC ラベル

最初の必要事項

- 2 機器の LAN インタフェースを以下に接続します。
 - クロスオーバー ネットワーク ケーブルを使用して PC のネットワーク カード (二点間)、または
 - 標準 LAN ケーブルを使用してハブまたはスイッチ



TCP/IP パラメータ コンフィグレーション

TCP/IP パラメータ コンフィグレーション

ネットワーク環境で正しく操作するには、有効な TCP/IP ネットワークパ ラメータを使用して LAN インターフェイスを設定する必要があります。こ のパラメータにはつぎのものが含まれます。

- IP アドレス
- サブネットマスク
- デフォルトゲートウェイ

以下の方法で TCP/IP パラメータを設定できます。

- ネットワークベースの BOOTP サーバからパラメータを自動的に要求する (いわゆる Bootstrap プロトコルを使用)。
- ネットワークベースの DHCP サーバからパラメータを自動的に要求する (いわゆる動的ホストコンフィグレーションプロトコルを使用)。この 方法にはオンボード LAN モジュールまたは G1369C LAN インターフェイ スカードが必要です(『「セットアップ(DHCP)」64ページ』を参照)。
- Telnet を使用して、手動でパラメータを設定する。

LAN インターフェイスは、複数の初期化モードを区別します。初期化モード(短縮形「init mode」)により、電源投入後に TCP/IP パラメータを有効にする方法が定義されます。パラメータは Bootp サイクルもしくは非揮発性メモリから派生、または既知のデフォルト値によって初期化されます。初期化モードは、コンフィグレーションスイッチで選択します。『58 ページ表 11』を参照してください。

コンフィグレーション スイッチ

コンフィグレーション スイッチ

機器背面左側でコンフィグレーション スイッチにアクセスできます。

Agilent 1220 Infinity LC は、スイッチ 7 と 8 の設定が **ON** の状態で出 荷されます。これは、機器がデフォルトの固定 IP アドレスに設定されて いることを意味します。192.168.254.11

注記 LAN を設定するには、SW1 および SW2 を OFF に設定する必要があります。

表 10 出荷時のデフォルト設定

初期化(「Init」)モード	デフォルトを用いて、スイッチ 7 と 8 の 設定を On に切り替えます。
リンクコンフィグレーション	自動ネゴシエーションで決定される スピードおよびデュプレックスモード

注記

G4294B では、機器の LAN アクセスを設定するためには、DAD メイン ボード上 のコンフィグレーション スイッチを使用する必要があります。1220 Infinity LC メイン ボードはすべてオフに設定されている必要があります。 4 LAN コンフィグレーション 初期化モード選択

初期化モード選択

以下の初期化(init)モードを選択できます。

表 11 初期化モード切り替え

	SW 6	SW 7	SW 8	初期化モード
ON	オフ	オフ	オフ	Bootp
	オフ	オフ	オン	Bootp および保存
	オフ	オン	オフ	保存されたパラメータを使用
	オフ	オン	オン	デフォルトを使用
	オン	オフ	オフ	DHCP ¹

¹ ファームウェア B.06.40 以降が必要。オンボード LAN なしのモジュールについては、G1369C LAN イン ターフェイス カードを参照

Bootp

初期化モード Bootp が選択された場合、モジュールは Bootp サーバから パラメータのダウンロードを試みます。取得されたパラメータは、すぐに 有効なパラメータになります。モジュールの非揮発性メモリには保存され ません。そのため、パラメータはモジュールの電源をオン / オフすると失 われます。



Bootp および保存

Bootp および保存が選択された場合、Bootp サーバから取得されたパラ メータはすぐに有効なパラメータになります。さらに、モジュールの非揮 発性メモリに保存されます。そのため、電源を一旦切って入れ直した後も 利用可能です。これにより、モジュールの Bootp once コンフィグレー ションの一種が有効になります。

例:ネットワーク内で常に Bootp サーバがアクティブであることをユー ザーが望まない場合があります。しかし一方で、Bootp 以外のコンフィグ レーションメソッドを持っていない場合があります。この場合、Bootp サーバを一時的に起動して、初期化モードの Bootp および保存を使用して モジュールの電源を入れ、Bootp サイクルが完了するのを待ち、Bootp サーバを終了して、モジュールの電源を切ります。次に、初期化モードの 保存されたパラメータを使用を選択して、モジュールの電源を再度入れま す。今後は、その 1 回の Bootp サイクルで取得されたパラメータを使用 して、モジュールに対する TCP/IP 接続を確立できます。



非揮発性メモリへの書き込みは時間を要するため、初期化モードの Bootp お よび保存の使用には注意が必要です。そのため、電源を入れるごとにモジュー ルが Bootp サーバからパラメータを取得するようにする場合、初期化モード は Bootp が推奨されます。

注記



保存されたパラメータを使用

初期化モードの保存されたパラメータを使用が選択された場合、パラメー タはモジュールの非揮発性メモリから取得されます。これらのパラメータ を使用して、TCP/IP 接続が確立されます。パラメータは、説明したメソッ ドの1 つで事前に設定されています。



図 13 保存されたパラメータを使用(原則)

デフォルトを使用

デフォルトを使用が選択された場合、工場出荷時のパラメータが取得され ます。これらのパラメータにより、追加設定を行うことなく LAN インター フェイスへの TCP/IP 接続が有効になります。『61 ページ 表 12』を参照し てください。



注記

LAN のデフォルトアドレスを使用するとネットワークに問題が生じる恐れがあります。注意して、すぐに有効なアドレスに変更してください。

初期化モード選択

IP アドレス:	192. 168. 254. 11
サブネットマスク:	255. 255. 255. 0
デフォルトゲートウェイ	指定なし

表 12 デフォルトパラメータを使用

デフォルト IP アドレスは、いわゆるローカルアドレスのため、ネット ワーク機器で経路指定されません。そのため、PC とモジュールは同じサブ ネット内に存在する必要があります。

ユーザーはデフォルト IP アドレスを使用して Telnet セッションを開き、 モジュールの非揮発性メモリに保存されたパラメータを変更できます。そ の後セッションを閉じ、初期化モードに保存されたパラメータを使用を選 択して電源を再び入れ、新しいパラメータを使用して TCP/IP 接続を確立 できます。

モジュールが、LAN から分離されて、PC に直接配線されている場合 (クロ スオーバーケーブルまたはローカルハブなどを使用)、ユーザーはデフォ ルトパラメータを簡単に保存して TCP/IP 接続を確立できます。

注記

デフォルトを使用モードでは、モジュールのメモリーに保存されたパラメータ は自動的には消去されません。ユーザーが変更しなければ、保存されたパラ メータを使用モードに切り替えた後も、それらをまだ使用できます。

動的ホストコンフィグレーションプロトコル(DHCP)

動的ホストコンフィグレーションプロトコル(DHCP)

一般情報 (DHCP)

動的ホストコンフィグレーションプロトコル (DHCP) は IP ネットワーク 上で使用される自動コンフィグレーションプロトコルです。DHCP 機能は、 オンボード LAN インターフェイスおよび「B」ファームウェア (B. 06.40 以降)を搭載したすべての Agilent HPLC モジュールで利用できます。

- G1314D/E/F VWD
- G1315C/D DAD
- G1365C/D MWD
- G4212A/B DAD
- G4220A/B バイナリポンプ
- G1369C LAN インターフェイスカード
- 1120/1220 LC システム

初期化モード「DHCP」が選択された場合、カードは DHCP サーバからパラ メータのダウンロードを試みます。取得されたパラメータは、すぐに有効 なパラメータになります。カードの非揮発性メモリには保存されません。

ネットワークパラメータの要求に加えて、カードは DHCP サーバにホスト 名を送信します。ホスト名はカードの MAC アドレスと同じになっています (0030d3177321 など)。ホスト名 / アドレス情報は DHCP サーバによってド メイン名サーバに送信されます。カードはホスト名解決のためのサービス (NetBIOS など)を一切提供しません。



図 15 DHCP (原則)

動的ホストコンフィグレーションプロトコル (DHCP)

注記

- 1 DHCP サーバがホスト名情報で DNS サーバを更新するまでにはしばらく時間 がかかることがあります。
- **2** DNS のサフィックス (0030d3177321. country. company. com など)を使用してホスト名を完全に限定する必要がある場合があります。
- **3** DHCP サーバがカードの提示したホスト名を拒否し、以下のローカルの命名 規則に従った名前を割り当てる場合があります。

動的ホストコンフィグレーションプロトコル (DHCP)

セットアップ (DHCP)

- **必要なソフト** スタックのモジュールには少なくともセット A. 06. 34 のファームウェアおよ ウェア: び前述のモジュール B. 06. 40 以降が必要です(同一のファームウェアセット のものである必要があります)。
 - 1 LAN インターフェイスの MAC アドレス (G1369C LAN インターフェイス カードまたはメインボードに付属)をメモしてください。MAC アドレス はカードのラベルまたはメインボードの背面に記載されています (0030d3177321 など)。

インスタントパイロットの MAC アドレスは、LAN セクションの 詳細で 確認できます。

	System Info	
Property	Value	
Contro	oller : DE12345678 (G4208A)	Relead
Main Revision	B.02.12 [0001]	Reibau
DAI	D : DE64260019 (G1315D)	
Main Revision	B.06.41 [0002]	
Resident Revison	B.06.40 [0007]	Print
On-time	3d 01:33h	
Installed Options	Dhcp	
LAN TCP/IP Mode	DHCP	
LAN TCP/IP Address	130.168.132.219	
LAN MAC Address	0030D314F89E	
Board ID	TYPE=G1315-66565, SER=MAC, REV=AC, MFG=	
Lamp	2140-0820 : 848728	
Cell	no info	
	Ţ	Exit
Information on each m	odule.	10:08

図 16 インスタントバイロットの LAN 設定

動的ホストコンフィグレーションプロトコル(DHCP)

- 2 前述のモジュールの G1369C LAN インターフェイスカードおよびメイン ボードのコンフィグレーションスイッチを DHCP にします。
- 表 13 G1369C LAN インターフェイスカード (カードのコンフィグレー ションスイッチ)

SW 4	SW 5	SW 6	SW 7	SW 8	初期化モード
オン	オフ	オフ	オフ	オフ	DHCP

表 14 1120/1220 を含む LC モジュール (機器背面のコンフィグレー ションスイッチ)

SW 6	SW 7	SW 8	初期化モード
オン	オフ	オフ	DHCP

- **3** LAN インターフェイスのホストとなるモジュールの電源を入れます。
- 4 コントロールソフトウェア (Agilent ChemStation、LabAdvisor、ファームウェア更新ツールなど)を設定し、MAC アドレス(例: 0030d3177321)をホスト名に使用します。

コントロールソフトウェアに LC システムが表示されます(『「一般情報 (DHCP)」62ページ』の注を参照)。

リンクコンフィグレーション選択

リンクコンフィグレーション選択

LAN インターフェイスは、フルまたはハーフデュプレックスモードで 10 または 100 Mbps の動作をサポートしています。多くの場合で、ネット ワークスイッチまたはハブなどのネットワーク機器を接続する場合にフル デュプレックスがサポートされ、IEEE 802.3u 自動ネゴシエーション仕様 をサポートします。

自動ネゴシエーションをサポートしないネットワーク機器を接続する場合、 LAN インターフェイスはそれ自体を 10 または 100 Mbps ハーフデュプ レックス動作に設定します。

たとえば、非ネゴシエーション 10 Mbps ハブに接続それる場合、LAN イン ターフェイスは 10 Mbps ハーフデュプレックスで動作するように自動的に 設定されます。

モジュールが自動ネゴシエーションを介してネットワークに接続できない 場合、モジュールのリンクコンフィグレーションスイッチを手動で設定で きます。

表 15 リンクコンフィグレーションのスイッチ

	SW 3	SW 4	SW 5	リンク コンフィグレーション
	オフ	_	-	自動ネゴシエーションで決定される スピードおよびデュプレックスモー ド
1 2 3 4 5 6 7 8	オン	オフ	オフ	10 Mbps、ハーフデュプレックスに マニュアル設定
	オン	オフ	オン	10 Mbps、フルデュプレックスにマ ニュアル設定
	オン	オン	オフ	100 Mbps、ハーフデュプレックスに マニュアル設定
	オン	オン	オン	100 Mbps、フルデュプレックスにマ ニュアル設定

BootP を使用した自動コンフィグレーション

BootP を使用した自動コンフィグレーション

注記	この章で示したすべての例が、ご使用の環境で動作するとは限りません。自身の IP アドレス、サブネットマスクアドレス、ゲートウェイアドレスが必要です。
注記	検出器のコンフィグレーションスイッチを確実に正しく設定するようにしてく ださい。設定は、BootP または BootP および保存 のいずれかにする必要があ ります (『58 ページ 表 11』を参照)。
注記	ネットワークに接続された検出器の電源がオフになっていることを確認してく ださい。
注記	Agilent BootP サービスプログラムがご使用の PC にインストールされていな い場合、Agilent ChemStation DVD にある BootP フォルダからインストールし てください。

Agilent BootP サービスについて

Agilent BootP サービスを使用して、LAN インターフェイスに IP アドレ スを割り当てます。

Agilent BootP サービスは ChemStation DVD で提供されます。Agilent BootP サービスは、LAN のサーバまたは PC にインストールされ、LAN 上 のアジレント機器の IP アドレスを一元管理します。BootP サービスは TCP/IP ネットワークプロトコルを実行している必要があり、DHCP サーバ を実行することはできません。

BootP を使用した自動コンフィグレーション

BootP サービスの使用

機器の電源を入れると、機器の LAN インターフェイスが IP アドレスまた はホスト名の要求を送信し、そのハードウェア MAC アドレスを識別子とし て提供します。Agilent BootP サービスはこの要求に応答して、直前に定 義された IP アドレスと、ハードウェア MAC アドレスに関連付けられたホ スト名を要求元の機器に渡します。

機器はその IP アドレスとホスト名を受け取り、電源が入っている限りこ の IP アドレスを維持します。機器の電源を切ると IP アドレスが失われ るため、機器に電源を入れるたびに Agilent BootP サービスを実行しなけ ればなりません。Agilent BootP サービスをバックグラウンドで実行して いる場合は、機器は電源投入時にその IP アドレスを受け取ります。

Agilent LAN インターフェイスは IP アドレスを保存するように設定でき るため、電源を入れ直しても IP アドレスは失われません。

状況:LAN 通信を確立できない場合

BootP サービスとの LAN 通信が確立できない場合は、PC で次の内容を確認します。

- BootP サービスが起動しているかどうか。BootP のインストール時は、 サービスは自動的に起動されません。
- ファイアウォールが BootP サービスをブロックしているかどうか。 BootP サービスを例外として追加します。
- LAN インターフェイスが「保存またはデフォルト」モードではなく BootP モードを使用しているかどうか。

BootP を使用した自動コンフィグレーション

BootP サービスのインストール

Agilent BootP サービスをインストールし、構成する前に、必ずコン ピュータと機器の IP アドレスを手元に用意してください。

- 1 管理者として、または管理者権限を持つ他のユーザーとしてログオンします。
- **2** すべての Windows プログラムを閉じます。
- 3 Agilent ChemStation ソフトウェア DVD をドライブに挿入します。セットアッププログラムが自動的に起動したら、キャンセル をクリックして停止します。
- 4 Windows Explorer を開きます。
- **5** Agilent ChemStation DVD の BootP ディレクトリに移動し、 BootPPackage.msi をダブルクリックします。
- 6 必要に応じて、タスクバーの Agilent BootP Service... アイコンをク リックします。
- 7 Agilent BootP Service Setup Wizard の Welcome 画面が表示されます。 次へ をクリックします。
- End-User License Agreement 画面が表示されます。条件を読んだ後、承認して 次へ をクリックします。
- 9 Destination Folder 選択画面が表示されます。BootP をデフォルトの フォルダにインストールするか、または Browse をクリックして別の場 所を選択します。次へ をクリックします。

デフォルトのインストール場所は次のとおりです。

C:\Program Files\Agilent\BootPService\

10 Install をクリックしてインストールを開始します。

BootP を使用した自動コンフィグレーション

11 ファイルがロードされます。完了すると、BootP 設定 画面が表示されま す。

ootP Tab File: :\Documents and Settings\ Create Tab File	All Users\Application Data\4	gilent\BootP\Ta	hFile	
NDocuments and Settings:	All Users\Application Data\4	gilent\BootP\Ta	hEile	
Create Tab File	1		01110	
	Edit BootP Address	es		
Logging				
🗖 Do you want to log boo	otP requests?			
BootP Log File:				
C:\Documents and Setting	IsVAII UsersVApplication Data	\Agilent\BootP\L	.ogFile	
Default Settings				
Subnot mode	0 0 0 0	-		
Subnet mask.				
Gateway:	0.0.0.0	-		
		OK	Cancel	Help

図 17 [BootP 設定] 画面

12 スクリーンの初期設定 部分では、もしわかっているなら、サブネット マスクとゲートウェイを入力できます。

デフォルトが使用できます:

- デフォルト サブネットマスクは、255.255.255.0 です。
- デフォルトゲートウェイは、192.168.254.11 です。
- **13** BootP 設定 画面で OK をクリックします。Agilent BootP Service Setup 画面に操作の完了が表示されます。
- **14** Finish をクリックし、Agilent BootP Service Setup 画面を終了しま す。
- 15 DVD をドライブから取り外します。

これでインストールは完了です。

16 Windows ® サービスの BootP サービスを開始します: Windows ® デスクトップで、コンピュータアイコンをクリックし、管理 > サービスとアプリケーション > サービスを選択します。Select the Agilent BootP サービス を選択して、開始をクリックします。

BootP を使用した自動コンフィグレーション

MAC アドレスを決定する 2 つの方法

ロギングを有効にし、BootP を使用して MAC アドレスを検出する

MAC アドレスを表示するには、Do you want to log BootP requests? チェックボックスをオンにします。

- 1 [BootP 設定] を、スタート > すべてのプログラム > Agilent BootP Service > EditBootPSettings の順に選択して開きます。
- **2** BootP 設定... で Do you want to log BootP requests? をオンにし、 ロギングを有効にします。

[- Logging
	✓ Do you want to log bootP requests? BootP Log File:
	C:\Documents and Settings\All Users\Application Data\Agilent\BootP\LogFile

図 18 BootP ロギングの有効化

ログファイルは次の場所にあります。

C:¥Documents and Settings¥All Users¥Application Data¥Agilent¥ BootP¥LogFile

ここには、BootP から構成情報を要求する各デバイスの MAC アドレスエ ントリが含まれます。

- 3 OK をクリックして値を保存するか、キャンセル をクリックして破棄し ます。編集が終了します。
- 4 BootP の設定(EditBootPSettings など)を変更するたびに、BootP サービスを停止または開始して変更を承認する必要があります。 『「Agilent BootP サービスの停止」76ページ』または『「Agilent BootP サービスの再起動」77ページ』を参照してください。
- 5 機器の構成後に Do you want to log BootP requests? ボックスをオフ にします。オフにしないと、ディスクスペースはすぐにログファイルで いっぱいになります。

BootP を使用した自動コンフィグレーション

LAN インタフェース カードラベルから直接 MAC アドレスを決定

- 1 機器の電源を切ります。
- 2 ラベルから MAC アドレスを読み込んで記録します。

MAC アドレスはモジュール背面のラベルに印刷されています。バーコードの下のコロン(:)の右側に記載されている、通常は AD から始まる番号が MAC アドレスです。『54ページ 図 9』および 『55ページ 図 10』を参照してください。

3 機器の電源を入れます。
BootP を使用した自動コンフィグレーション

Agilent BootP サービスによる IP アドレスの割り当て

Agilent BootP サービスでは、機器のハードウェア MAC アドレスを IP ア ドレスに割り当てます。

BootP サービスを使用した機器の MAC アドレスの確認

- 1 機器の電源を一旦切り、入れ直します。
- 2 機器の自己診断が終了したら、メモ帳を使用して BootP サービスのログ ファイルを開きます。
 - ログファイルのデフォルトの場所は C:¥Documents and Settings¥All Users¥Application Data¥Agilent¥BootP¥LogFile です。
 - ログファイルは、開いていると更新されません。

内容は次のようなものになります。

02/25/10 15:30:49 PM

Status:BootP Request received at outermost layer

Status:BootP Request received from hardware address:0010835675AC

Error:Hardware address not found in BootPTAB:0010835675AC

Status:BootP Request finished processing at outermost layer

- **3** ハードウェア (MAC) アドレス (0010835675AC など) を記録します。
- **4** Error は、MAC アドレスが IP アドレスに割り当てられておらず、タブ ファイルにこのエントリがないことを示します。IP アドレスが割り当 てられると、MAC アドレスはこのタブファイルに保存されます。
- 5 別の機器の電源を入れる前にログファイルを閉じます。
- 6 機器の構成後に Do you want to log BootP requests? ボックスをオフ にして、ログファイルで過度のディスクスペースを使用しないようにし ます。

BootP を使用した自動コンフィグレーション

BootP を使用して各機器をネットワークに追加

- 1 スタート > すべてのプログラム > Agilent BootP Service の順に選択 し、Edit BootP Settings を選択します。[BootP 設定] 画面が表示さ れます。
- 2 すべての機器が追加されたら、Do you want to log BootP requests? を オフにします。

機器の構成が完了したら **Do you want to log BootP requests?** ボック スをオフにします。オフにしないと、ディスクスペースはすぐにログ ファイルでいっぱいになります。

- 3 Edit BootP Addresses... を選択します。Edit BootP Addresses 画面が 表示されます。
- 4 追加... をクリックします。Add BootP Entry 画面が表示されます。

Mac Address	
Host Name	
IP Address	· · ·
Comment	
Subnet Mask	255 . 255 . 255 . 0
Gateway	· · ·
<u>0</u> K	<u>Cancel</u> <u>H</u> elp

図 19 BootP ロギングの有効化

BootP を使用した自動コンフィグレーション

- 5 機器について次の項目を入力します。
 - ・ MAC アドレス
 - ホスト名。希望のホスト名を入力します。
 ホスト名の先頭は「アルファベット」でなければなりません(GC6890 など)。
 - IP アドレス
 - コメント(オプション)
 - サブネットマスク
 - ゲートウェイアドレス(オプション)

入力した構成情報はタブファイルに保存されます。

- 6 OK をクリックします。
- 7 Close を押して Edit BootP Addresses を終了します。
- 8 OK をクリックして BootP 設定 を終了します。
- BootPの設定([EditBootPSettings] など)を変更するたびに、BootP サービスを停止または開始して変更を承認する必要があります。
 『「Agilent BootP サービスの停止」76ページ』または『「Agilent BootP サービスの再起動」77ページ』を参照してください。
- 10機器の電源を一旦切り、入れ直します。

または

IP アドレスを変更した場合は、機器の電源を一旦切って入れ直すこと により変更を有効にします。

11 PING ユーテリティを使用して、接続性を確認するため、コマンドウィ ンドウを開いて、以下を打ち込みます:

例えば、Ping 192.168.254.11。

タブファイルは、次の場所に位置します。

C:¥Documents and Settings¥All Users¥Application Data¥Agilent¥ BootP¥TabFile

BootP を使用した自動コンフィグレーション

Agilent BootP サービスによる機器の IP アドレスの変更

PC を再起動すると、Agilent BootP サービスは自動的に起動します。 Agilent BootP サービスの設定を変更するには、サービスを停止し、変更 を行い、サービスを再起動する必要があります。

Agilent BootP サービスの停止

1 Windows のコントロールパネルで、管理ツール > サービス の順に選択 します。サービス 画面が表示されます。



- 図 20 Windows の [サービス] 画面
- 2 Agilent BootP Service を右クリックします。
- 3 停止 を選択します。
- 4 サービスおよび管理ツール 画面を閉じます。

EditBootPSettings での IP アドレスおよびその他のパラメータの 編集

- 1 スタート > すべてのプログラム > Agilent BootP Service の順に選択 し、Edit BootP Settings を選択します。BootP 設定 画面が表示され ます。
- **2** BootP 設定 画面を初めて開くと、インストール時のデフォルト設定が 表示されます。

BootP を使用した自動コンフィグレーション

3 Edit BootP Addresses... を押してタブファイルを編集します。

Hardware Address	Host Name	IP Address	Comment	Subnet Mask	Gateway
006000111999 005000222888	AgilentLC1 AgilentLC2	10.1.1.101 101.1.1.102	Agilent LC1 right Agilent LC2 left	255.255.255.0 255.255.255.0	0.0.0 0.0.0.0
q					

図 21 [Edit BootP Addresses] 画面

4 Edit BootP Addresses... 画面で、追加... をクリックして新しいエント リを作成するか、またはテーブルから既存の行を選択し、Modify... ま たは Delete を押して、IP アドレス、コメント、サブネットマスクなど をタブファイルで変更します。

IP アドレスを変更した場合は、機器の電源を一旦切って入れ直すことに より変更を有効にする必要があります。

- 5 Close を押して Edit BootP Addresses... を終了します。
- 6 [OK] をクリックして [BootP 設定] を終了します。

Agilent BootP サービスの再起動

- **1** Windows のコントロールパネルで、管理ツール > サービス の順に選択 します。サービス 画面が表示されます(『76ページ 図 20』を参照)。
- 2 Agilent BootP Service を右クリックし、開始 を選択します。
- 3 サービスおよび管理ツール 画面を閉じます。

Bootp を使用した設定の永久保存

Bootp を使用した設定の永久保存

Bootp を使用してモジュールのパラメータを変更する場合、以下の指示に 従ってください。

- 1 モジュールの電源を切ります。
- 2 モジュールの設定スイッチの設定を「Bootp および保存」モードに変更 します(『58ページ表11』を参照)。
- **3** Agilent Bootp サービスを起動して、そのウィンドウを開きます。
- 4 必要に応じて、既存のコンフィグレーションを使用してモジュールのパ ラメータを修正します。
- **5** [OK] を押して、[Bootp マネージャ] を終了します。
- 6 ここでモジュールの電源を入れ、[Bootp サーバ] ウィンドウを表示させ ます。しばらくした後、Agilent Bootp サービスにより LAN インタ フェースからのリクエストが表示されます。パラメータは、ここでモ ジュールの非揮発性メモリに恒久的に保存されます。
- **7** Agilent Bootp サービスを閉じて、モジュールの電源を切ります。
- 8 モジュールの設定スイッチの設定を「保存されたパラメータを使用」モー ドに変更します(『58ページ表11』を参照)。
- 9 モジュールの電源をオン / オフします。これで、Agilent Bootp サービス 無しで LAN 経由でアクセスできます。

LAN コンフィグレーション 4 マニュアルコンフィグレーション

マニュアルコンフィグレーション

マニュアルコンフィグレーションでは、モジュールの非揮発性メモリに保存されたパラメータだけが変更されます。現在のアクティブパラメータに影響を及ぼすことはありません。そのため、いつでもマニュアルコンフィグレーションを実行できます。保存されたパラメータをアクティブパラメータにするには電源のオン / オフが必須で、これにより初期化モード選択スイッチが有効となります。



4 LAN コンフィグレーション マニュアルコンフィグレーション

Telnet の使用

モジュールへの TCP/IP 接続が可能な場合(いずれかの方法で TCP/IP パ ラメータ設定されている場合)はいつでも、Telnet セッションを開くこと でパラメータを変更できます。

- 1 Windows の [スタート] ボタンをクリックして、[名前を指定して実行] を選択して、システム (DOS) プロンプトウィンドウを開きます。「cmd」 と打ち込み、[OK] をクリックします。
- 2 システム (DOS) プロンプトで以下のように入力します。
 - c:¥>telnet <IP アドレス>または
 - c:¥>telnet < ホスト名 >



図 23 Telnet - セッションの開始

ここで〈IP address〉は、Bootp サイクル、ハンドヘルド コントローラ を使用したコンフィグレーションセッション、あるいはデフォルトの IP アドレスから割り当てられたアドレスになります(『「コンフィグ レーション スイッチ」57ページ』を参照)。

接続が確立された場合、モジュールは以下のように応答します。



図 24 モジュールとの接続が確立されます。

3?と打ち込み、Enterを押すと、使用可能なコマンドを確認できます。

C:\WINDOWS\system32\cmd.exe - telnet 134.40.27.95		
Agilent Technologies G ≻?	L315C PP00000024	
command syntax	description	
? / ip <x.x.x.x> sm <x.x.x.x> gw <x.x.x.x> exit ></x.x.x.x></x.x.x.x></x.x.x.x>	display help info display current LAN settings set IP Address set Subnet Mask set Default Gateway exit shell	

図 25 Telnet コマンド

表 16 Telnet コマンド

値	説明
?	構文およびコマンドの説明を表示します
/	現在の LAN 設定を表示します
ip <x. x="" x.=""></x.>	新しい IP アドレスを設定します
sm <x. x="" x.=""></x.>	新しいサブネットマスクを設定します
gw <x. x="" x.=""></x.>	新しいデフォルトゲートウェイを設定します
exit	シェルを終了して、すべての変更を保存します

4 パラメータを変更するには、以下のスタイルに従ってください。

パラメータ値、たとえば:

ip 134.40.27.230

次に [Enter] を押します。ここでパラメータは定義しているコンフィグ レーションパラメータを参照して、値はパラメータに割り当ててある定 義を参照します。各パラメータ入力後、改行します。

マニュアルコンフィグレーション

5「/」を使用して、[Enter]を押すと、現在の設定が一覧表示されます。

C:\WINDOWS\s	ystem32\cmd.exe - telnet 134.40.27.95
>∕ LAN Status Page	8
MAC Address	: 0030D30A0838
Init Mode	: Using Stored
TCP/IP Propert: - active - IP Address Subnet Mask Def. Gateway	ies : 134.40.27.95 : 255.255.248.0 : 134.40.24.1
TCP/IP Status	: Ready
Controllers >_	: no connections

LAN インターフェイス、MAC アドレス、 MAC アドレス、初期化モード 初期化モード - [保存されたパラメータ を使用] アクティブな TCP/IP 設定 TCP/IP ステータス - ここでは レディ コントロールソフトウェア (Agilent ChemStation など)を使用した PC との 接続 - ここでは接続されていません

- 図 26 Telnet 「保存されたパラメータを 使用」モードでの現在の設定
 - 6 IP アドレスを変更して (この例では、134.40.27.99)、「/」を打ち込み、 現在の設定を一覧表示します。

95

C:\WINDOWS\s	system32\cmd.exe - telnet 134.40.27
>ip 134.40.27.	.99
LAN Status Pag	ye
MAC Address	: 0030D30A0838
Init Mode	: Using Stored
TCP/IP Propert	ties
IP Address	: 134.40.27.95
Subnet Mask	: 255.255.248.0
Def. Gateway	: 134.40.24.1
- stored -	
IP Address	: 134.40.27.99
Subnet Mask	: 255.255.248.0
Def. Gateway	: 134.40.24.1
TCP/IP Status	: Ready
Controllers	: no connections
図 2/ leine	at - IP 設定の変更

IP 設定の変更 初期化モード - [保存されたパラメータ を使用] アクティブな TCP/IP 設定 非揮発性メモリに保存された TCP/IP 設定

コントロールソフトウェア (Agilent ChemStation など)を使用した PC との 接続 - ここでは接続されていません

マニュアルコンフィグレーション

コンフィグレーションパラメータを打ち込み終ったら、
 exit を打ち込み、Enter を押し、パラメータを保存しながら終了します。



図 28 Telnet セッションの終了

注記

ここで初期化モードスイッチが「保存されたパラメータを使用」モードに変更 されると、モジュールが再起動された際に装置は保存された設定を取得しま す。上記の例では、134.40.27.99 になります。 4 LAN コンフィグレーション マニュアルコンフィグレーション



1220 Infinity LC

5

溶媒送液システムの説明

概要 86 デガッサ 87 操作原理 88 圧縮率補正 92 可変ストローク容量 94 ポンプの使用 95

この章では、溶媒送液システム(ポンプとオプションのデガッサ) の操作原理の概要を説明します。



5 溶媒送液システムの説明 概要



ポンプは、2 チャンネルのデュアルプランジャ直列型設計に基づく製品で、 溶媒送液システムに要求されるすべての機能を装備しています。最大 600 bar の圧力を発生できる 1 つのポンプアセンブリによって、溶媒のメ タリングと高圧側への送液が実行されます。

デガッサ内で溶媒が脱気され、高速定量バルブの低圧側で溶媒を混合しま す。デュアル チャンネルグラジエント ポンプには、内蔵デュアルチャン ネルオンラインデガッサが含まれます。Agilent 1220 Infinity LC のアイ ソクラティック ポンプにはデガッサは搭載されていません。

ポンプアセンブリには、注入口バルブとアウトレットバルブの付いたポン プヘッドが含まれます。ダンピングユニットは、2 つの プランジャチャン バの間に接続されます。ポンプアウトレットは、ポンプヘッドのプライミ ングに便利なように、PTFE フリットを含むパージバルブが取り付けられて います。

デガッサ

デュアル チャンネル グラジエント ポンプには、内蔵オンラインデガッサ が付属しています。ポンプの電源を入れると、流量が 0 mL/min に設定さ れていても、デガッサの電源は自動的に入ります。2 つのチャンネルの真 空チャンバに、75 Torr (100 mbar) の一定した真空ができます。溶媒は、 PTFE AF チューブ中を、真空チャンバ内の内部容量 1.5 mL/ チャンネルで 流れます。

5 **溶媒送液システムの説明** 操作原理

操作原理

液体は、溶媒ボトルからデガッサを通って DCGV へ流れ、そこから注入口 バルブへ流れます。ポンプアセンブリは、2 つの実質的に同一のプラン ジャ / チャンバユニットから構成されています。プランジャ / チャンバの 両ユニットは、ボールスクリュードライブと往復運動するサファイアプラ ンジャ 1 本を搭載したポンプヘッドから構成されています。

サーボ制御方式の可変磁気抵抗モータが、2 つのボールスクリュードライ ブをそれぞれ逆方向に駆動します。2 つのボールスクリュードライブ用の ギアの円周はそれぞれ異なる(2:1 の比)ため、第 1 プランジャは、第 2 プランジャの 2 倍の速度で動きます。溶媒は、ポンプヘッドの最下部近く から入って、ポンプヘッドの最上部から出ます。プランジャの外径はポン プヘッドチャンバの内径より小さいため、溶媒がそのすき間を充填します。 第 1 プランジャのストローク量は、流量に応じて 20 - 100 µL で変化し ます。すべての流量は、マイクロプロセッサによって、1 µL/min ~ 10 mL/min の範囲でコントロールされます。第 1 プランジャ / チャンバユ ニットの注入口は、注入口バルブに接続され、その注入口バルブは、第 1 プランジャポンプユニットに溶媒を吸引できるように開閉されます。

第 1 プランジャ / チャンバユニットの出口は、アウトレットボールバルブ とダンピングユニットを通して、第 2 プランジャ / チャンバユニットの注 入口に接続されます。次に、パージバルブアセンブリの出口は、クロマト グラフシステムに接続されます。

溶媒送液システムの説明 5 操作原理



ポンプの基本原理 図 29

ポンプの電源を入れると、ポンプは初期化手順を実行して、第1プラン ジャの上死点を決定します。第1プランジャは、チャンバの機械的停止位 置までゆっくりと上向きに動き、そこからあらかじめ指定した距離だけ下 向きに動きます。コントローラは、このプランジャ位置をメモリに保存し ます。この初期化後、ポンプは設定パラメータを使用して動作を開始しま す。注入口バルブが開き、プランジャが下向きに動いて、第1 チャンバに 溶媒を吸引します。同時に、第2プランジャが上向きに動き、システムへ 送液します。コントローラが(流量に応じて)ストローク長を定義した後、 ドライブのモータが停止し、注入口バルブが閉じます。モータが逆方向に 回転して、保存された上限値に達するまで第1 プランジャを上向きに動か し、同時に第2プランジャを下向きに動かします。次に、一連の動作が再 び開始され、プランジャを上限値と下限値の間で上下に動かします。第1 プランジャが上向きに動いている間、チャンバ内の溶媒はアウトレット ボールバルブを介して、第2 チャンバに押し出されます。第2 プラン ジャは、第1プランジャによって押し出された容量の半分を吸引します。 残り半分の容量は、直接システム内に送液されます。第 1 プランジャの吸 引ストローク中に、第2プランジャは吸引した量をシステムに送液しま す。

溶媒ボトル A、B からの溶媒の混合比に応じて、コントローラは吸引スト ロークの長さを分割します。分割された各部分で、グラジエントバルブは 指定の溶媒チャンネルをポンプ入力に接続します。

表 17 アイソクラティックポンプの詳細

デッド ボリューム	800 - 1100 µL(背圧により異なる)
移動相と接触する部品	の材質
ポンプヘッド	SST、金、サファイア、セラミック
アクティブインレッ トバルブ	SST、金、サファイア、ルビー、セラミック、PTFE
アウトレットバルブ	SST、金、サファイア、ルビー
アダプタ	SST、金
パージバルブ	SST、金、PTFE、セラミック、PEEK
デガッサチャンバ	TFE/PDD 共重合体、FEP、PEEK、PPS

表	18	グラジエン	トポン	プの詳細
---	----	-------	-----	------

ディレイ ボリューム	800 - 1100 µL (背圧により異なる)
移動相と接触する部品の	D材質
MCGV	PTFE
ポンプヘッド	SST、金、サファイア、セラミック
アクティブインレッ トバルブ	SST、金、サファイア、ルビー、セラミック、PTFE
アウトレットバルブ	SST、金、サファイア、ルビー
アダプタ	SST、金
パージバルブ	SST、金、PTFE、セラミック、PEEK
ダンパーユニット	金、SST
デガッサチャンバ	TFE/PDD 共重合体、FEP、PEEK、PPS

圧縮率補正

圧縮率補正の原理

システム内の背圧が変化した場合(カラムの老化など)、使用する溶媒の圧 縮率の違いが、リテンションタイムの安定性に影響を与えます。この影響 を最小限に抑えるために、ポンプは溶媒のタイプに応じて流量安定性を最 適化する圧縮率補正機能を装備しています。圧縮率補正値はデフォルト値 に設定されており、ユーザーインタフェースを使用して変更できます。

圧縮率補正なしで、第1プランジャのストローク中に以下のことが生じま す。プランジャチャンバの圧力が上昇し、チャンバ内の容量が圧縮されま す。これは、背圧と溶媒タイプにより異なります。このため、圧縮された 容量の分だけ、システムに送液される容量が減少してしまいます。

圧縮率の値を設定しておけば、プロセッサが補正容量を自動的に計算しま す。補正容量は、システム内の背圧と選択した圧縮率によって決まります。 通常のストローク容量にこの補正容量を加算することによって、すでに説 明した第1プランジャの送液ストローク中に起こる容量の損失を補正しま す。

圧縮率補正設定の最適化

圧縮率補正のデフォルト設定は、46・10⁻⁶ /bar です。この設定値は平均値 です。通常の条件下では、このデフォルト設定で、圧力の脈動を大部分の アプリケーションやすべてのグラジェント分析に十分な値(システム圧力 の1%未満)に抑えます。高感度検出器を用いたアプリケーションでは、 さまざまな溶媒の値を用いることで圧縮率設定を最適化できます。また、 アイソクラティック混合液を使用する際に、溶媒が圧縮率の表に含まれて いない場合、そして初期設定がご使用のアプリケーションに対して十分で はない場合は、以下の手順で圧縮率設定を最適化できます。

注記 混合溶媒を使用する場合、その混合液に使用される純粋な溶媒の圧縮率値を補 間したり、その他の計算を適用しても、混合液の圧縮率を計算することは不可 能です。このような場合、以下の実験的手順を適用して、圧縮率設定を最適化 する必要があります。 1 必要な流量でポンプを起動します。

- 2 最適化手順を開始する前に、流量が安定している必要があります。脱気した溶媒だけを使用します。耐圧テストでシステムの気密を確認します。
- 3 ご使用のポンプを、圧力とリップル(%)をモニタリングできるコント ロールソフトウェアに接続する必要があります。
- 4 圧力リップルが最少になる圧縮率補正値が、使用する溶媒の組成に最適 な設定値です。

溶媒(原液)	压縮率(1.10 ⁻⁶ /bar)
アセトン	126
アセトニトリル	115
ベンゼン	95
四塩化炭素	110
クロロホルム	100
シクロヘキサン	118
エタノール	114
エチルアセテート	104
ヘプタン	120
ヘキサン	150
イソブタノール	100
イソプロパノール	100
メタノール	120
1- プロパノール	100
トルエン	87
水	46

表 19 溶媒圧縮率

5 溶媒送液システムの説明 可変ストローク容量

可変ストローク容量

ポンプチャンバ内の容量の圧縮によって、ポンプのプランジャストローク のたびに小さな圧力の脈動が発生し、ポンプの流量リップルに影響を与え ます。圧力の脈動の振幅は、主にストローク量と使用中の溶媒の圧縮率補 正値によって決まります。ストローク量が小さいと、同じ流速ではスト ローク量が大きい場合よりも、振幅の小さい圧力の脈動が生じます。さら に、圧力の脈動の周波数が高くなります。したがって、定量結果に対する 流量変動の影響が小さくなります。

グラジエントモードでは、ストローク量が小さいほど、流量リップルが小 さくなり、混合によるリップルが改善されます。

ポンプは、プロセッサ制御方式のスピンドルシステムを使用してプラン ジャを駆動します。通常のストローク量は、選択した流量に合わせて最適 化されます。流量が小さい場合は、ストローク量も小さくなり、流量が大 きい場合はストローク量も大きくなります。

ポンプのストローク量が AUTO モードに設定されている場合、ストローク は流量に合わせて最適化されます。ストローク量をこれより大きくするこ とも可能ですが、この変更はお勧めできません。

ポンプの使用

Agilent 1220 Infinity LC ポンプをうまく使用するための ヒント

- Agilent 1120 Infinity LC ポンプに塩溶液と有機溶媒を使用する際は、 塩溶液を下部グラジエントバルブポートの1 つに接続し、有機溶媒を上 部グラジエントバルブポートの1 つに接続することをお勧めします。有 機チャンネルを塩溶液チャンネルのすぐ上にするのがベストです。すべ ての DCGV チャンネルを水で定期的にフラッシングし、バルブポートの 塩堆積物をできるだけ取り除くことをお勧めします。
- ポンプを運転する前に、特に一定期間ポンプの電源が切られていた場合 (終夜など)、および揮発性溶媒混合液をチャンネルに使用した場合は、
 2 倍以上の量(3 mL)でデガッサ(オプション)をフラッシュします。
- 溶媒インレットフィルタの詰まりを防止します(溶媒インレットフィル タなしでポンプを決して使用しないでください)。藻の繁殖を防止する 必要があります。
- ・ 定期的にパージバルブフリットとカラムフリットを確認します。パージバルブフリットの詰まりは、その表面に黒または黄色の層ができていたり、圧力が10 bar 以上(パージバルブを開いて5 mL/min の流量で蒸留水を送液する場合)になることで特定できます。
- 低流量(たとえば、0.2 mL/min)でポンプを使用する場合、すべての 1/16 インチフィッティングに漏れの兆候がないかを確認します。
- ポンプシールを交換するときは、パージバルブフリットも交換してください。
- バッファ溶液を使用した場合、システムの電源を切る前に水でフラッシュします。
- プランジャのシールを交換する場合、ポンププランジャに傷がないか確認します。プランジャに傷があると、微量リークの原因になり、シールの寿命を縮める可能性があります。
- プランジャシールの交換後、馴らし作業に従ってシステムを加圧します。

溶媒フィルタの詰まり防止

溶媒が汚れていたり、溶媒ボトル内に藻が増殖したりすると、溶媒フィル タの寿命が短くなるばかりでなく、ポンプの性能に悪影響を与えます。水 性溶媒またはリン酸緩衝液(pH4 - 7)を使用する場合は、特に注意が必 要です。溶媒フィルタの寿命を延ばし、ポンプの性能を維持するために、 次の注意に従ってください。

- ・藻の繁殖を遅らせるために、可能であれば、茶色の溶媒ボトルを使用してください。
- 溶媒は、藻を除去するフィルタまたは被膜で濾過してください。
- ・ 溶媒は2 日ごとに交換するか、濾過し直してください。
- アプリケーションで可能であれば、溶媒に 0.0001 0.001 M のアジ化 ナトリウムを添加してください。
- 溶媒の上にアルゴン層を形成すると藻の発生が防げます。
- ・ 溶媒ボトルを直射日光にさらさないでください。

注記

溶媒フィルタを取り付けずにシステムを使用しないでください。



1220 Infinity LC

6

注入システムの説明

マニュアルインジェクタ 98
インジェクションシール 99
サンプルの注入 99
ニードル 101
オート サンプラ 102
サンプリング動作 103
注入動作 104
サンプリングユニット 106
ニードルドライブ 107
アナリティカルヘッド 107
注入バルブ 108
トランスポート アセンブリ 109
オート サンプラでサポートされるトレイ 110
バイアルとキャップの選択 111

この章では、注入システム(マニュアルインジェクタおよびオート サンプラ)の操作原理の概要を説明します。



マニュアルインジェクタ

Agilent 1220 Infinity LC マニュアルインジェクタでは、Rheodyne、6 ポートサンプルインジェクションバルブ (5067-4202)を使用しています。 サンプルはバルブ正面のインジェクションポートを通って外部の 20 µL の サンプルループに充填されます。バルブには PEEK[™] インジェクションシー ルが施されています。ステータには make-before-break passage が設けら れているため、バルブの位置が 注入ポジションからロード ポジションに 切り替えられた場合や、元の位置に戻された場合でも、溶媒の流れは中断 されません。





マニュアルインジェクタ

インジェクションシール

マニュアルインジェクタには、標準で PEEK[™] インジェクションシールが付 属しています。

サンプルの注入

警告

移動相の排出

100 µL 以上のサンプルループを使用する場合、サンプルループ中の 移動相が減圧するとニードルポートから移動相を排出できます。

→ 試薬メーカーから提供されている取扱説明書および安全データ シートの記載に従って、適切な安全手順(ゴーグル、安全手袋、 防護衣など)を守ってください。特に、有毒溶媒や有害な溶媒を 使用する場合は、注意してください。

ロードポジション

ロードポジション(『100ページ 図 31』を参照)で、ポンプを直接カラム (ポート 2 と 3 が接続された)に接続し、ニードルポートをサンプル ループに接続します。精度を高めるためには、2 ~ 3 のサンプルループ 容量(さらに高い精度が必要な場合はさらに多く)をニードルポートから 注入する必要があります。サンプルでループを満たし、余分なサンプルは ポート 6 に接続されたベントチューブから放出されます。

6 注入システムの説明

マニュアルインジェクタ



図 31 ロードポジション

注入ポジション

注入ポジション(『100ページ 図 32』を参照)で、ポンプをサンプルルー プ(ポート1 ~ 2 が接続された)に接続します。すべてのサンプルを ループからカラムへ洗い出します。ニードルポートをベントチューブ (ポート5)に接続します。





外径 0.028 インチ (22 ゲージ) x 長さ 2 インチ、電気的なテーパ加工無 し、90 °の先端様式 (正方形チップ)のニードルを使用します。

オート サンプラ

オート サンプラには 3 種類のサイズのサンプルラックを利用できます。 標準フルサイズのラックは、100 × 1.8 mL バイアルを収容し、一方 2 個のハーフサイズラックのスペースは、40 × 1.8 mL および 15 × 6 mL バイアルをそれぞれ収容します。オート サンプラには同時に 2 枚のハーフサイズラックトレイを設置できます。アナリティカルヘッド デバイスにより、下限 0.1 - 100 µL の量から注入できます。

オート サンプラの移送機構は、X-Z-シータ運動を使用して、バイアルの ピックアップと返却を最適化します。バイアルはグリッパアームでピック アップされ、サンプリング ユニット下の位置に置かれます。グリッパの移 送機構とサンプリングユニットはモーター駆動です。正しく動作するよう に、動作は光学センサと光学エンコーダでモニタリングされます。キャ リーオーバを最小限に抑えるために、注入後にメタリングデバイスは必ず フラッシュされます。

6 ポートの注入バルブユニット(使用するのは 5 ポートのみ)は、高速 ハイブリッドステッパモータ駆動です。サンプリング動作中は、バルブユ ニットはオート サンプラをバイパスして、ポンプからカラムへフローを直 接接続します。注入および分析時は、バルブユニットはオート サンプラに フローを向けて、サンプルが完全にカラム内に注入されるようにし、その 後、次のサンプリング動作が始まる前にメタリングユニットとニードルか らサンプル残留物を除去するようにします。

サンプリング動作

サンプリングシーケンス中のオートサンプラコンポーネントの動作は、プ ロセッサによって常時モニタリングされています。このプロセッサによっ て、各動作のタイムウィンドウと物理的な移動範囲が定義されています。 サンプリング動作の特定ステップを正常に終了できないと、エラーメッ セージが生成されます。

サンプリング動作中、インジェクションバルブによって溶媒はオートサン プラからバイパスされます。グリッパアームで固定サンプルラックからサ ンプルバイアルを選択します。グリッパアームにより、注入ニードル下に サンプルバイアルを置きます。メタリングデバイスで必要なサンプル量を サンプルループに吸引します。サンプリング動作の終了時点で注入バルブ がメインパスポジションに戻ると、サンプルがカラムに注入されます。

サンプリング動作は、次の順序で実行されます。

- 1 インジェクションバルブがバイパスポジションに切り換えられます。
- 2 メタリングデバイスのプランジャが初期化ポジションに移動します。
- 3 グリッパアームでバイアルを選択します。同時に、ニードルはシートの 外に上昇します。
- 4 グリッパアームにより、ニードル下にバイアルを置きます。
- 5 ニードルがバイアルの中に下降します。
- 6 メタリングデバイスにより、設定されたサンプル量を吸引します。
- 7 ニードルがバイアルから上昇します。
- 8 自動ニードル洗浄が選択されると、グリッパアームによりサンプルバイ アルを取り替え、ニードル下に洗浄バイアルを置き、バイアルの中に ニードルを下げ、その後洗浄バイアルの外にニードルを上昇させます。
- 9 安全フラップが所定の位置にあるか、グリッパアームで確認します。
- 10 グリッパアームでバイアルを交換します。同時に、ニードルはシートの 中に下がります。
- 11 インジェクションバルブはメインパス位置に切り替わります。

注入動作

注入動作の開始前、および分析中は、インジェクションバルブはメインパ スポジションです。インジェクションバルブがこのポジションにあると、 移動相はオートサンプラのメタリングデバイス、サンプルループ、および ニードル内を送液されます。これにより、サンプルに触れた部分がすべて 測定中にフラッシュされ、キャリーオーバを最小限に抑えます。



図 33 メインパスポジション

サンプリング動作開始時、バルブユニットはバイパスポジションに切り替わります。ポンプから送られた移動相は、ポート1のバルブユニットに入り、ポート6を通ってカラムに直接流れます。



図 34 バイパスポジション

次にニードルを上昇させ、ニードルの下にバイアルを置きます。ニードル はバイアルの中に下がり、メタリングユニットによりサンプルをサンプル ループに吸引します。



メタリングユニットにより必要量のサンプルをサンプルループに吸引する と、ニードルを上昇し、バイアルをサンプルトレイに戻します。ニードル はニードルシートの中に下がり、インジェクションバルブはメインパスポ ジションに戻り、カラムにサンプルを流します。



サンプリングユニット

サンプリングユニットは、ニードルドライブ、メタリングドライブ、イン ジェクションバルブの3つのメインアセンブリから構成されています。

注記

交換用サンプリングユニットには、インジェクションバルブとメタリングヘッ ドアセンブリは含まれません。

ニードルドライブ

ニードルの動作は、ギアベルトでスピンドルアセンブリに接続されたス テッパモータにより駆動されます。モータの回転運動は、スピンドルアセ ンブリのドライブユニットで直線運動に変換されます。ニードルの上端位 置と下端位置はサンプリングユニットフレックスボードの反射センサで検 知されますが、ニードルのバイアル内ポジションはニードルセンサの上端 ポジションからモータステップを計数することで測定します。

アナリティカルヘッド

アナリティカルヘッドは、ギアベルトでドライブのシャフトに接続された ステッパモータによって駆動します。スピンドルの回転運動は、スピンド ル上の駆動ナットによって、直線運動に変換されます。サファイアのプラ ンジャは、駆動ナットによってスプリングに押し付けられ、アナリティカ ルヘッドに入ります。プランジャの底部は駆動ナットの大きなベアリング 上にあるため、プランジャは常に中心の位置にあります。プランジャは、 セラミックリングによってアナリティカルヘッドに導かれます。プラン ジャのホームポジションは、サンプリング ユニットのフレックスボード上 にある赤外線センサによって検出されます。一方、サンプル量は、ホーム ポジションからのステップ数を計数することで決定します。プランジャが 後ろに下がる(スプリングによって駆動される)ことによって、サンプル がバイアルから吸引されます。

	標準 (100 µL)
ステップ数	15000
分解能	7 nL/モータのステップ
最大ストローク	100 µ L
圧力限界値	600 bar
プランジャの材質	サファイア

表 20 アナリティカルヘッドの技術データ

注入バルブ

2 ポジション 6 ポート注入バルブは、1 つのステッパモータによって駆動 します。使用されるのは、6 つのポートのうち、5 つだけです(ポート 3 は使用されません)。ステッパモータの動きは、レバー/スライダ機構に よって注入バルブに伝達されます。注入バルブの切り換え(バイパスポジ ションとメインパスポジション)は、2 つのマイクロスイッチによってモ ニタリングされます。

内部のコンポーネントを交換した後のバルブの調整は不要です。

	標準
モーターのタイプ	4 V、1.2 A ステッパモータ
シールの材質	Vespel™(Tefzel™使用可能)
ポート数	6
スイッチング時間	$< 150 {\rm \ ms}$

表 21 注入バルブの技術データ
トランスポート アセンブリ

トランスポートユニットは、X 軸スライド(左右動作)、Z 軸アーム(上下 動作)、グリッパアセンブリ(回転とバイアルのグリップ)から構成されて います。



図 37 トランスポートアセンブリ

トランスポートアセンブリでは、サンプル バイアル移送用のグリッパアセ ンブリを正確に位置決めするために、閉ループモードで駆動する 4 つのス テッパモータを使用します。モーターの回転運動は、ドライブスピンドル に接続されたギアベルトにより直線運動(X 軸と Z 軸)に変換されます。 グリッパアセンブリの回転(シータ軸)は、ギアベルトと一連のギアによ りモーターから伝えられます。グリッパフィンガーの開閉は、グリッパア センブリ内部の遊星ギアにギアベルトでリンクされたステッパモータによ り駆動されます。 ステッパモータの位置は、ステッパモータのハウジングに設置された光学 エンコーダで決定されます。エンコーダにより、モーターの位置を継続し てモニタリングし、ポジションエラーを自動的に修正します(たとえば、 バイアルトレイ内にバイアルを収納する際に、グリッパが偶発的にズレた 場合)。可動コンポーネントの初期化位置は、フレックスボードに設置され た反射センサで検知されます。これらの位置は、プロセッサが実際のモー ター位置を計算するために使用されます。さらに、トレイ認識用の 6 つの 反射センサがアセンブリ前面のフレックスボードに設置されています。

オート サンプラでサポートされるトレイ

部品番号	説明
G1313-44510	トレイ、100 x2 mL バイアル
G1313-44513	ハーフトレイ、15 x6 mL バイアル
G1313-44512	ハーフトレイ、40 x2 mL バイアル

ハーフ トレイの組み合わせ

ハーフ トレイは任意のタイプを組み合わせた取り付けが可能で、2 mL- お よび 6 mL- バイアルを同時に使用できます。

バイアルポジションの番号付け

標準の 100 - バイアルトレイには、1 ~ 100 のバイアルポジションがあ ります。しかし、2 枚のハーフ トレイを使用する場合は、番号付けの規 則が少し異なります。右側のハーフトレイのバイアルポジションは以下に 示すようにポジション 101 から始まります。

左側 40 - ポジションのトレイ:1-40

- 左側 15 ポジションのトレイ:1 15
- 右側 40 ポジションのトレイ: 101 140
- 右側 15 ポジションのトレイ: 101 115

バイアルとキャップの選択

信頼性の高い操作を行うために、Agilent 1220 Infinity LC オート サン プラで使用されるバイアルは、テーパー形の肩やバイアル本体より幅広い キャップではないようにする必要があります。下表に部品番号と共に示し たバイアルとキャップは、Agilent 1220 Infinity LC オート サンプラを 用いて最小 15000 回注入したテストに合格しました。

クリンプバイアル

部品番号	説明
5181-3375	クリンプバイアル、2 mL、透明ガラス、100 個入
5183-4491	クリンプバイアル、2 mL、透明ガラス、1000 個入
5182-0543	クリンプバイアル、2 mL、透明ガラス、ラベル付き、 100 個入
5183-4492	クリンプバイアル、2 mL、透明ガラス、ラベル付き、 1000 個入
5183-4494	クリンプバイアル、2 mL、透明ガラス、ラベル付き、 100 個入(シラン処理)
5181-3376	クリンプバイアル、2 mL、褐色ガラス、ラベル付き、 100 個入
5183-4493	クリンプバイアル、2 mL、褐色ガラス、ラベル付き、 1000 個入
5183-4495	クリンプバイアル、2 mL、褐色ガラス、ラベル付き、 100 個入(シラン処理)
5182-0567	クリンプバイアル、1 mL、ポリプロピレン製、広口、 100 個入
5183-4496	クリンプバイアル、1 mL、ポリプロピレン製、広口、 1000 個入(シラン処理)
9301-0978	クリンプバイアル、0.3 mL、ポリプロピレン製、広口、 1000 個入

6 注入システムの説明 オート サンプラ

スナップバイアル

部品番号	説明
5182-0544	スナップバイアル、2 mL、透明ガラス、100 個入
5183-4504	スナップバイアル、2 mL、透明ガラス、1000 個入
5183-4507	スナップバイアル、2 mL、透明ガラス、100 個入(シラン 処理)
5182-0546	スナップバイアル、2 mL、透明ガラス、ラベル付き、 100 個入
5183-4505	スナップバイアル、2 mL、透明ガラス、ラベル付き、 1000 個入
5183-4508	スナップバイアル、2 mL、透明ガラス、ラベル付き、 100 個入(シラン処理)
5182-0545	スナップバイアル、2 mL、褐色ガラス、ラベル付き、 100 個入
5183-4506	スナップバイアル、2 mL、褐色ガラス、ラベル付き、 1000 個入
5183-4509	スナップバイアル、2 mL、褐色ガラス、ラベル付き、 100 個入(シラン処理)

スクリューバイアル

部品番号	説明
5182-0714	スクリューキャップバイアル、2 mL、透明ガラス、 100 個入
5183-2067	スクリューバイアル、2 mL、透明ガラス、1000 個入
5183-2070	スクリューバイアル、2 mL、透明ガラス、100 個入 (シラン処理)
5182-0715	スクリューバイアル、2 mL、透明ガラス、ラベル付き、 100 個入
5183-2068	スクリューバイアル、2 mL、透明ガラス、ラベル付き、 1000 個入
5183-2071	スクリューバイアル、2 mL、透明ガラス、ラベル付き、 100 個入(シラン処理)
5182-0716	スクリューキャップバイアル、2 mL、褐色ガラス、ラベル 付き、100 個入
5183-2069	スクリューバイアル、2 mL、褐色ガラス、ラベル付き、 1000 個入
5183-2072	スクリューバイアル、2 mL、褐色ガラス、ラベル付き、 100 個入(シラン処理)

6 注入システムの説明

オート サンプラ

クリンプキャップ

部品番号 説明

- 5181-1210 クリンプキャップ、銀色アルミ、セプタム (透明 PTFE/赤色ラバー)、100 個入
- 5183-4498クリンプキャップ、銀色アルミ、セプタム
(透明 PTFE/ 赤色ラバー)、1000 個入
- 5181-1215 クリンプキャップ、青色アルミ、セプタム (透明 PTFE/赤色ラバー)、100 個入
- 5181-1216クリンプキャップ、緑色アルミ、セプタム
(透明 PTFE/赤色ラバー)、100 個入
- 5181-1217 クリンプキャップ、赤色アルミ、セプタム (透明 PTFE/赤色ラバー)、100 個入

スナップキャップ

部品番号 説明

- 5182-0550スナップキャップ、透明ポリプロピレン、セプタム
(透明 PTFE/赤色ラバー)、100 個入
- 5182-3458スナップキャップ、青色ポリプロピレン、セプタム
(透明 PTFE/赤色ラバー)、100 個入
- 5182-3457スナップキャップ、緑色ポリプロピレン、セプタム
(透明 PTFE/赤色ラバー)、100 個入
- 5182-3459スナップキャップ、赤色ポリプロピレン、セプタム(透明 PTFE/赤色ラバー)、100 個入

スクリューキャップ

部品番号 説明

- 5182-0717スクリューキャップ、青色ポリプロピレン、セプタム
(透明 PTFE/赤色ラバー)、100 個入
- 5182-0718スクリューキャップ、緑色ポリプロピレン、セプタム
(透明 PTFE/赤色ラバー)、100 個入
- 5182-0719スクリューキャップ、赤色ポリプロピレン、セプタム
(透明 PTFE/赤色ラバー)、100 個入
- 5182-0720スクリューキャップ、青色ポリプロピレン、セプタム
(透明 PTFE/ シリコン)、100 個入
- 5182-0721スクリューキャップ、緑色ポリプロピレン、セプタム
(透明 PTFE/シリコン)、100 個入
- 5182-0722スクリューキャップ、赤色ポリプロピレン、セプタム
(透明 PTFE/シリコン)、100 個入

6 注入システムの説明 オート サンプラ



この章では、カラムオーブンの操作原理の概要を説明します。



7 カラムオーブンの説明 カラムオーブン

カラムオーブン

カラムオーブンは、2 つの温度センサを搭載した抵抗ヒーターマットに基づき、カラム領域全体の温度を一定にします。内蔵の過熱温度遮断ヒューズにより、オーバーヒートを抑制します。

オーブンキャピラリの内部容量は 6 µL です。

カラムの長さの最大は、25 cm (10 inch) です。

操作範囲は、室温の 5 ° 以上、少なくとも 10 ° C から最高 60 ° C ま で、指定流速は最大 5 mL/min (60 ° C) です。

注記 フロント カバーが必ず閉じた状態で正しいカラム温度になるように、フロン ト カバーを開けた状態でカラムオーブンを操作しないでください。オーブン 分離の片方は、フロント カバーの内側に固定されています。



1220 Infinity LC

8 検出器の説明

検出器タイプ 120 Agilent 1220 Infinity LC 可変波長検出器 (VWD) 121 検出器 121 Agilent 1220 Infinity ダイオードアレイ検出器 (DAD) 122 検出器の概要 122 光学系 123 ピーク幅(レスポンスタイム) 126 サンプルおよびリファレンス波長および帯域幅 128 スリット幅 131 スペクトル取り込みの最適化(DAD のみ) 133 ネガティブ吸収のマージン 133 選択性の最適化 134 スペクトル設定(DAD のみ) 137 適切なカラムとフローセルの組合せ 140

この章では、検出器の操作原理の概要を説明します。





検出器タイプ

Agilent 1220 Infinity LC システムには、2 つの異なる検出器タイプを 使用できます。

- 可変波長検出器 (VWD、G4286B、G4288B/C、G4290B/C で使用)、G1314F
 VWD の光学ユニット
- ダイオードアレイ検出器 (DAD、G4294B で使用)、G1315C DAD の光学ユ ニット

Agilent 1220 Infinity LC 可変波長検出器 (VWD)

Agilent 1220 Infinity LC 可変波長検出器 (VWD)

検出器

Agilent 1220 Infinity LC 可変波長型検出器は、優れた光学的性能を発揮 し、GLP に準拠し、メンテナンスが容易に行えるように設計されています。

- 重水素ランプを内蔵し、波長範囲 190 600 nm で最大強度と最低検出 限界を実現、
- オプションのフローセルカートリッジ(標準:10 mm14 µL、高圧: 10 mm14 µL、マイクロ:3 mm2 µL、セミマイクロ:6 mm5 µL)が使用可 能で、アプリケーションの必要に応じて使用できます、
- ランプとフローセルは前面から容易にアクセスでき、すばやく交換でき、
- 内蔵の酸化ホルミウム フィルタにより、波長真度の確認をすばやく行 えます。

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)

検出器の概要

本検出器は、優れた光学的性能を発揮し、GLP に準拠し、メンテナンスが 容易に行えるように設計されています。本検出器には、以下のような特徴 があります。

- 80 Hz (超) 高速 LC アプリケーション用のデータ取込速度
- すべてのフローセルおよび UV- ランプに RFID (無線認識) タグがあり、 アセンブリの情報の追跡が可能です。
- RFID (無線認識) タグの付いた長寿命の重水素およびタングステン ランプにより、最大強度と最低検出限界を実現できる波長範囲は 190 950 nm
- 同時に最大 8 波長に対し、感度喪失なし
- 下限 1 16 nm のプログラム式スリットで、感度、直線性、スペクトル 分解の完全な最適化
- RFID (無線認識) のあるオプションのフローセルカートリッジ (標準 10 mm13 µL、セミマイクロ 6 mm5 µL、マイクロ 3 mm2 µL、80 nL、 500 nL、10 mm、高圧 10 mm1.7 µL およびプレップーセル) が、アプリ ケーションの必要に応じて使用可能です。
- ランプとフローセルは前面から容易にアクセスでき、すばやく交換可能
- 内蔵の酸化ホルミウム フィルタにより、波長真度をすばやく確認
- ・ 内蔵温度コントロール (ETC) により、ベースラインの安定性を向上
- ・ 温度およびランプ電圧のモニタリング用追加診断シグナル

仕様については、『「Agilent 1220 Infinity LC DAD の性能仕様」27ページ』を参照してください。

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)

光学系

検出器の光学系は、下図に示す通りです。光源は、紫外線(UV) 波長範囲 用の重水素アーク放電ランプと、可視光線(VIS) および短波近赤外 (SWNIR) 範囲用のタングステン ランプの組み合わせです。タングステン ランプのフィラメントのイメージは、特別なバックアクセス ランプ設計に より、重水素ランプの放電開口部に焦点が合います。これにより、両光源 が光学的に組み合わされ、ソース レンズに対し軸を共有できます。アクロ マート(ソース レンズ)は、フローセルを通って単一の集中光線を形成し ます。各セル ルームとランプはクォーツ窓で仕切られており、クォーツ窓 はクリーンアップまたは交換できます。分光器では、光りはホログラ フィー グレーティングによりダイオードアレイに分散します。これによっ て、すべての波長情報に同時にアクセスすることが可能になります。



図 38 検出器の光学系

ランプ: UV- 波長範囲の光源は、シャインスルー型開口部をもつ重水素ランプです。 低圧重水素ガス内でのプラズマ放電によって、ランプは波長 190 nm から 約 800 nm までの範囲の光を放出します。可視光線と SWNIR の波長範囲用 光源は、低ノイズのタングステン ランプです。このランプは、470 ~ 950 nm の波長範囲で光を放出します。

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)

- **アクロマート** アクロマートは、両方のランプから光を受けると、光を収束させ、その (ソース ビームがフローセルを通過します。 レンズ)
- 酸化ホルミウム 酸化ホルミウム フィルタは、電気機械的に作動します。ホルミウムフィル フィルタ タ テスト中、フィルタは光路に移動します。
 - **セルサポート** セルサポートウィンドウのアセンブリは、ホルミウム フィルタエリアとフ **ウィンドウ** ローセルエリアを分離します。
- フローセルの コンパート メント
 光学ユニットにはフローセルのコンパートメントがあり、フローセルへの アクセスが容易です。オプションの種々のフローセルは、同一の簡単でシ ンプルな取り付けシステムを使用し、挿入することができます。フローセ ルは取り外して、フローセルに影響されずに検出器の光学的および電子的 パフォーマンスをチェックできます。
 - 分光器 分光器の材質はセラミック製で、温度効果を最小限に抑えます。分光器は、 分光器用レンズ、可変入射スリット、フロントエンド電子工学のグレー ティングおよびフォトダイオードアレイで構成されます。分光器用レンズ は、光ビームがフローセルを通過した後、光ビームを再収束させます。ダ イオードアレイのデータ間隔は、190 ~ 950 nm の波長範囲で、< 1 nm です。波長により、ナノメータ当り 1.0 から 1.25 ダイオードまで変動し ます (たとえば、0.8 から 1 nm ごとに 1 ダイオード)。

波長範囲が小さい場合は、小さな非線形は無視できます。波長範囲が 190 ~ 950 nm の場合は、すべての範囲で波長真度を達成するためには新しい アプローチが必要です。各分光器は個別にキャリブレーションされます。 キャリブレーション データは、分光器の EEPROM に保存されます。これら のデータに基づき、内蔵プロセッサが、データポイント間は直線間隔 (1.0、2.0、…) で、吸光度データを計算します。その結果、優れた波長真 度と機器間の再現性が実現します。

- **可変入射** マイクロスリットシステムは、シリコンの機械的特性にバルク微細加工技 スリット 術の精密構造を組み合わせて開発されました。必要な光学的機能 – スリッ
- **システム** トおよびシャッターはシンプルでコンパクトな構成要素を組み合わせてい ます。スリット幅は、機器のマイクロプロセッサによって直接制御され、 メソッドパラメータとして設定することができます。
- **グレーティング** 凹型ホログラフィー グレーティングを使用して、分散とスペクトル画像解 析を達成します。グレーティングは、光ビームを各波長に分光し、その光 をフォトダイオードアレイ上に反射します。

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)

ダイオード アレイ アレイ ア上に組み込んだ制御回路から構成されています。波長範囲 190 ~ 950 nm では、データ間隔は< 1 nm です。

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)

ピーク幅(レスポンスタイム)

レスポンスタイムとは、フローセル内の吸光度の突然の変化に検出器のシ グナルがいかに速く追随するかを示すものです。検出器は、デジタルフィ ルタを使用して、レスポンスタイムをクロマトグラムのピーク幅に適合さ せます。フィルタは、ピーク面積やピーク対称度には影響しません。正し く設定されていれば、フィルタによりベースラインノイズは大幅に減少し ます(『126ページ図39』)が、ピーク高さもわずかに減少します。さら に、フィルタによりデータ速度を遅くすることで、ピークの積分と表示を 最適化し、クロマトグラムとスペクトルを保存するのに必要なディスクス ペースを小さくすることができます。



図 39 レスポンスタイムのシグナルとノイズへの影響

『127 ページ 表 22』に、検出器のフィルタの選択肢を一覧表示します。最 適な結果を得るには、ピーク幅を、クロマトグラム中の目的ピークにでき るだけ近くなるように設定します。レスポンスタイムは、ピーク幅の約 1/3 に設定します。結果として、ピーク高さは 5 % 未満減少し、ピーク分 散が 5 % 未満増加します。検出器のピーク幅設定を小さくすると、ピーク 高さが 5 % 未満増加しますが、レスポンスタイムが 1/2 になるとベース ラインノイズが 1.4 倍に増加します。ピーク幅(レスポンスタイム)を 推奨設定の 2 倍にすると(オーバーフィルタリング)、ピーク高さは約 20 % 低くなり、ベースラインノイズは 1.4 分の 1 になります。シグナ ル/ノイズ比はベストになりますが、ピーク分解能に影響が出ます。

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)

ピーク幅[分]	レスポンスタイム [秒]	データ速度 [Hz]
< 0.0025	0.025	80
> 0.0025	0.05	80
> 0.005	0.1	40
> 0.01	0.2	20
> 0.03	0.5	10
> 0.05	1.0	5
> 0.10	2.0	2.5
> 0.20	4.0	1.25
> 0.40	8.0	0.62
> 0.85	16.0	0.31

表 22 ピーク幅 - レスポンス タイム - データ速度

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)

サンプルおよびリファレンス波長および帯域幅

本検出器は、190 ~ 950 nm の波長で吸光度を同時に測定します。2 個の ランプにより、全波長範囲にわたって高感度です。重水素放電ランプが、 UV 範囲 (190 ~ 400 nm) 用のエネルギーを提供し、タングステン ランプ が、可視光線と短波近赤外用に 400 ~ 950 nm の光を放出します。

サンプル中の検体が未知の場合は、両方のランプを使用し、全波長範囲に わたってすべてのスペクトルを保存します。膨大なディスク空き容量が必 要となりますが、すべての情報を取込むことができます。スペクトルは、 ピークの純度チェックと定性に使用できます。また、スペクトル情報は、 クロマトグラフ シグナルの波長設定を最適化するためにも有用です。

本検出器は、以下のプロパティを基にして、分析時間に最大 8 シグナルを 取込み保存することができます。

- サンプル波長、サンプル帯域幅(BW)の波長バンドの中心、およびオプション
- リファレンス波長、リファレンス帯域幅の波長バンドの中心。

シグナルは、サンプル波長バンドの平均吸光度からリファレンス波長バンドの平均吸光度を差し引いて、時間に対するの一連のデータポイントで構成されます。

検出器デフォルトメソッドのシグナル A は、サンプル 250、100、リファ レンス 360、100 に設定されています。すなわち、200 ~ 300 nm の平均吸 光度から、300 ~ 400 nm の平均吸光度を差し引いたものです。すべての 検体は、200 ~ 300 nm の範囲で、300 ~ 400 nm の範囲における吸光度よ り高い吸光度を示すので、このシグナルで、UV 吸光のある多くの化合物を 検出できます。

多くの化合物がそのスペクトラムで吸光帯を示します。『129 ページ 図 40』は、アニス酸のスペクトルの例です。

アニス酸の検出下限濃度で最適化するには、サンプル波長を吸光帯のピーク(252 nm)に設定し、サンプル帯域幅を吸光帯の幅(30 nm)に設定します。360、100のリファレンスは適切です。アニス酸は、この範囲で吸収を持ちません。

高濃度で分析する場合は、サンプル波長をスペクトルの谷に設定すること により 1.5 AU 以上での良好な直線性が得られます。アニス酸の場合は 225 nm です。

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)



図 40 波長設定の最適化

帯域幅を広くすると、広い波長範囲にわたって平均化されることにより、 ノイズが減少するという利点が得られます。帯域幅 4 nm に比べて、ベー スラインノイズは約 2.5 分の 1 になり、一方シグナルは帯域幅 4 nm の 場合の約 75 % になります。この例の場合、帯域幅 30 nm の場合の シグナ ル/ノイズ比は、帯域幅 4 nm の場合の 2 倍になります。



1220 Infinity LC

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)

検出器は、各波長について計算される吸光度値を平均化するので、広い帯 域幅を使用しても、直線性に悪影響を与えません。

ベースラインドリフト、および室温変動やグラジエント中の屈折率の変化 によるうねりを減少させるために、リファレンス波長の設定を強くお勧め します。

『130 ページ 図 42』に、PTH アミノ酸分析のベースライン ドリフト減少の 例を示します。リファレンス波長を使用しない場合、クロマトグラムはグ ラジエントによる屈折率の変化のために下方にドリフトします。リファレ ンス波長を使用することにより、このドリフトはほとんど完全に除去され ます。このテクニックを使用して、グラジエント分析でも低ピコモル範囲 で PTH-アミノ酸の定量を行うことができます。



図 42 PTH アミノ酸のグラジエント分析 (各 1 pmol)、リファレンス波長 を使用した場合と使用しない場合

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)

スリット幅

検出器には、分光器の入射部に可変スリットがあります。これは、種々の 分析内容に対応して、検出器を最適化させるのに有効な装備です。

狭いスリットを使用すると、吸光度スペクトルが微細構造を持っている検体で、高いスペクトル分解能が得られます。そのようなスペクトルの例として、ベンゼンの例を示します。5 つの主吸光帯(フィンガ)は、幅 2.5 nm で互いに 6 nm しか離れていません。



図 43 1、4、16 nm のスリット幅の場合のベンゼン(原理)

広いスリットの場合は、多くの光がフローセルを通過します。『132 ページ 図 44』に示すように、これによりベースラインノイズは低くなります。

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)



図 44 スリット幅のベースラインノイズに対する影響

ただし、スリット幅が広くなると、分光器の光学分解能(波長の違いを区別する能力)が減少します。フォトダイオードは、スリット幅によって決まる波長範囲内の光を受けます。スリット幅16 nm を使用するとベンゼンの微細なスペクトル構造が消失するのはこのためです。

一方化合物のスペクトルの急なスロープ部分の波長では、吸光度は濃度に 対して厳密な直線性を示しません。

ベンゼンのように微細構造を持ち吸光度が大きく変化する物質は、非常に まれです。

ほとんど場合、スペクトルの吸光帯の幅は、むしろアニス酸の場合のよう に 30 nm です(『129ページ 図 40』を参照してください)。

多くの場合、スリット幅 4 nm でベストな結果が得られます。

微細なスペクトル構造を持つ化合物を同定する場合、またはスペクトルの スロープ部分の波長で高濃度(> 1000 mAU)の定量を行う場合は、狭いス リット(1 または 2 nm)を使用します。一方、シグナルの帯域幅を広げる ことで、ベースラインノイズを小さくすることができます。(デジタル) 帯域幅を使用して吸光度を平均化するので、直線性に影響はありません。

サンプルが非常に低濃度である場合は、広い(8 または 16 nm)スリット を使用します。必ずスリット幅と同じ帯域幅のシグナルを使用してください。

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)

スペクトル取り込みの最適化(DAD のみ)

すべてのスペクトルを保存するには、大量のディスク空き容量が必要です。 メソッドの最適化時または未知サンプルを分析するときには、すべてのス ペクトルを取り込むことは有用です。しかし、既知のサンプルを多数分析 するときには、すべてのスペクトルを取り込むと、必要性の低いデータで 多くのデータ容量を取られてしまいます。本検出器では、データ量を減ら し、しかも必要なスペクトル情報は保持しておく機能があります。

スペクトルオプションについては、『138 ページ 表 23』を参照してください。

範囲

サンプル中の化合物が吸収する波長範囲に、純度チェックやライブラリ サーチに有益な情報が含まれます。スペクトル保存の波長範囲を減らすこ とによっても、ディスク空き容量を節約できます。

ステップ

大部分の物質は、広い吸光帯を持っています。吸光帯の幅あたり 5 ~ 10 のデータポイントでスペクトルが取込されている場合に、スペクトル表示、 ピーク純度とライブラリサーチの実施がベストになります。アニス酸(前 に使用した例)の場合は、4 nm のステップで十分です。ただし、ステップ を 2 nm にすると、スペクトルをより最適に表示します。

スレッショルド

ピーク検出器を設定します。ピークコントロール保存モードが選択されて いる場合は、スレッショルドより高いピークのスペクトルのみを保存しま す。

ネガティブ吸収のマージン

ベースラインがやや負側に(約-100 mAU)ドリフトする場合、バランス時 に検出器のゲイン調整を行います。特殊な場合、例えば、吸収を持つ溶媒 によるグラジエント分析を行う場合は、ベースラインがさらに負の値にド リフトすることがあります。

そのような場合に限り、ネガティブ吸収についてのマージンを増加させ、 A/D コンバータのオーバーフローを防止します。

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)

選択性の最適化

ピーク抑制による不分離ピークの定量化

クロマトグラフでは、しばしば 2 つの化合物が分離できずに溶出します。 従来の 2 波長検出器では、各々一方の化合物のみが吸収する波長がある場 合にのみ、各波長で互いに独立して化合物を検出し定量することができま す。ただし、これはごく少数の化合物にしか当てはまりません。

ダイオードアレイ技術によるデュアルチャネル検出器の場合は、両方の化 合物の吸収波長範囲が重なっている場合でも、2 つの化合物の定量化が可 能です。この手順は、ピーク抑制またはシグナル減算と呼ばれます。例と して、カフェインが存在するサンプルのヒドロクロロチアジドの分析につ いて説明します。生物学的サンプル中のヒドロクロロチアジドを分析する 場合は、必ずカフェインが存在するリスクがあり、ヒドロクロロチアジド のクロマトグラフに干渉する可能性があります。『134 ページ 図 45』のス ペクトルに示すように、ヒドロクロロチアジドは 222 nm においてベスト な感度に検出されますが、そこではカフェインも大きな吸光度を示します。 そのため、従来の可変波長検出器では、カフェインが存在する場合にはヒ ドロクロロチアジドを定量的に検出することは不可能です。



図 45 ピークサプレッションの場合の波長選択

ダイオードアレイに基づく UV-Vis 検出器を使用し、リファレンス波長を 正しく選択すれば、定量検出が可能です。カフェインを差し引くには、リ

1220 Infinity LC

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)

ファレンス波長を 282 nm に設定する必要があります。この波長で、222 nm におけるカフェインの吸光度と等しくなります。それぞれの吸光度値を 減算すれば、カフェイン吸光度が除去されます。同様にして、カフェイン を定量する場合はヒドロクロロチアジドを差し引くことができます。この 場合は、波長を 204 nm に設定し、リファレンス波長を 260 nm に設定し ます。『135 ページ 図 46』に、ピーク抑制テクニックによるクロマトグラ フの結果を示します。

ただし、その結果としては、感度が減少します。サンプルシグナルは、リファレンス波長における吸光度の比率だけシグナル波長の吸光度が減少します。感度は 10 ~ 30% 低くなります。



図 46 リファレンス波長を使用するピークサプレッション

比率クォリファイアによる化合物クラスの選択検出

比率クォリファイアを使用して、複雑なサンプル中のある特定のクラスの みを分析できます。たとえば、生物学的サンプル中の元の薬物とその代謝 物を分析するような場合です。もう 1 つの例は、プレカラムまたはポスト カラム誘導体化後の誘導体物質の選択的分析です。そのサンプル群に適し たシグナル比を指定することで、目的ピークのみを選択的にプロットする ことができます。比がユーザー指定比範囲を外れている間は、シグナル出 力はゼロのままです。比が範囲内に入ると、シグナル出力は標準の吸光度 に対応し、フラットなベースライン上に明らかなシングルピークとして出 力されます。『136 ページ 図 47』と 『136 ページ 図 48』に例を示しま す。

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)



図 47 レシオクォリファイアの波長選択



図 48 レシオクォリファイアによる選択性

4 成分の混合物において、ビフェニルのみが記録されています。その他の 3 つのピークは、比率クォリファイアの基準を満たさないために出力はゼ ロとなり、非表示になっています。特徴的な波長 249 nm (1₁) および 224 nm (1₂) が、『136 ページ 図 47』に示されているスペクトルから検出 されています。比率範囲は 2 ~ 2.4 (2.2 ± 10%) に設定されています。 249 nm と 224 nm の間の比率がこの範囲内である場合にのみ、シグナルが プロットされます。4 つのピークのうち、3 番目のピークのみが基準を満 たします (『136 ページ 図 48』)。他のピークはプロットされません。 Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)

スペクトル設定(DAD のみ)

スペクトル設定を変更するために開きます。

- 1 スペクトル設定を変更するには、[検出器設定シグナル]を選択します。
- 2 [スペクトル] セクションで、ドロップダウンリストをクリックして、 パラメータを選択します。『138 ページ 表 23』には可能なパラメータが 表示されます。
- 3 必要に応じて、範囲、ステップ幅、スレッショルドを変更します。

UE Set up DAD Signals	DAD Signals : System-2	DAD Signals : System-2		
🔓 Control	Signals	_ Signals		
📲 Configuration	Store Sample,Bw Reference,Bw	Store Sample,Bw Reference,Bw		
🔐 Run Recovery	<u>A</u> : 🗹 254 4 360 100 🗧 nm	<u>A</u> : 🗹 254 4 360 100 븆 nm		
	B: □ 254 16 360 100 븆 nm	B: □ 254 16 360 100 븆 nm		
ﷺ Not Ready Information 쉽 Help	C: □ 210 8 360 100 븆 nm	C: 🔲 210 8 360 100 🗦 nm		
	D: 🗖 230 16 360 100 🖨 nm	D: 🗖 230 16 360 100 🖨 nm		
	E: 🗖 280 16 360 100 불 nm	E: 🗖 280 16 360 100 븆 nm		
	F: 280 16 360 100 🕈 nm	F: 🔲 280 16 360 100 茾 nm		
	G: 🗖 280 16 360 100 🕈 nm	G: 🗌 280 16 360 100 불 nm		
	H: 🗖 280 16 360 100 븆 nm	H: 🗌 280 16 360 100 🕇 nm		
	Spectrum	- Spectrum		
	Store: Every 2nd spectrum 💌	Store: Every 2nd spectrum		
	None Range: Apex+Baselines Apex+Slopes+Baselines	Range: 190 to 400 nm		
	Step: All in peak Every 2nd spectrum	Step: 2.0 nm		
	Threshold:	Threshold: 1.00 mAU		

図 49 スペクトル設定

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)

表 23 スペクトル設定

保存	「シグナル A」のどのポイントでスペクトルが取得され 保存されるかを定義します。シグナル A は「ピークコ ントロールスペクトル採取」のコントロールに使用さ れ、他のシグナルはスペクトル取得に影響を及ぼしませ ん。
なし	スペクトルの取り込みはありません。
頂点 + ベース ライン	ピークの頂点とベースラインのスペクトルを取り込みま す。
頂点 + スロープ + ベースライン	ピークの頂点、ベースライン、アップスロープ、ダウン スロープのスペクトルを取り込みます。
ピーク内すべて	ピーク内のすべてのスペクトルを取り込みます。
注	上記の3 つのスペクトル取り込みタイプメソッドは、 ピークコントロールスペクトル取り込みとも呼ばれます DAD に設定したスレッショルドとピーク幅のパラメータ により、検出器ファームウェアがピークを検出します。 ピークコントロールのスペクトル保存を使用する場合 は、これらのパラメータを設定し、必要なピークをすべ て認識させるようにします。積分アルゴリズムにはま た、積分イベントで設定したスレッショルドとびピーク 幅パラメータによるピークの検出が含まれます。
2 スペクトル毎	[すべて] では、スペクトルは連続して取込されます が、1 つおきのスペクトルだけが保存され、他のスペク トルは削除されます。これにより、保存するデータ量を 減らすことができます。
すべて	スペクトルは、ピーク幅の設定に従い連続して取り込ま れます。ピーク幅あたり 8 つのスペクトルを取り込み ます。1 つのスペクトルの取込時間は、ピーク幅を 8 で割った値より少し小さい、つまり 0.01 秒以上から 2.55 秒以下となります。
注	シグナル A にピークがない場合には、スペクトルはあ りません。他のシグナルでピークが検出されていてもス ペクトルは取り込まれません。

Agilent 1220 Infinity ダイオードアレイ検出器 (DAD)

表 23 スペクトル設定

範囲	範囲は、スペクトル保存における波長範囲を定義しま す。 リミット値:上限値と下限値はいずれも 190 ~ 950 nm (1 nm 刻み)。上限値は下限値より少なくとも 2 nm 長波 長側でなければなりません。
ステップ	ステップは、スペクトル保存における波長の分解能を定 義します。 リミット値:0.10 ~ 100.00 nm (0.1 nm 刻み)。
スレッショルド	スレッショルドは、期待される最も小さいピーク高さ (mAU 単位) です。ピーク検出器では、検出感度の数値 より低いすべてのピークを無視し、スペクトルの保存は 行いません。 リミット値: 0.001 ~ 1000.00 mAU、0.001 mAU 刻み。 頂点 + ベースライン、頂点 + スロープ + ベースライ ン、ピーク内すべてのモードに使用可能

適切なカラムとフローセルの組合せ

適切なカラムとフローセルの組合せ

『140ページ 図 50』は、使用するカラムに適した推奨フローセルを示して います。複数の選択肢がある場合、大きいフローセルを使用するとベスト な検出限界が得られます。小さいフローセルを使用するとベストなピーク 分解能が得られます。

WWD 用のフローセルの選択

カラム長	標準的な ピーク幅	并	授フローセル		
<= 5 cm	0.025 min				
10 cm	0.05 min		セミマイクロ フローセル		
20 cm	0.1 min			標準 フローセル	
>= 40 cm	0.2 min				
	標準的な 流量	0.2 ml/min	$0.2 \sim 0.4 \ { m ml/n}$	inO.4 \sim 0.8 ml/m	inl \sim 5 ml/min
カラム内径		1.0 mm	2.1 mm	3.0 mm	4.6 mm

図 50 フローセルの選択

適切なカラムとフローセルの組合せ

標準的なカラム 長さ	標準的な ピーク幅	推奨フローセル							
$T \leq 5 \text{ cm}$	0.025 min	ミクロま; セミナノ	たは						
10 cm	0.05 min			セミミクロ フローセル	1				 高耐圧 フローセル、
20 cm	0.1 min					標準 フロー	・セル		圧力 100 bar 以上用
\geq 40 cm	0.2 min								•
	標準的な 流量	0.01 mL/min	0.2	0.20. mL/min	4	0.4 mL/mir	. 0.4	1 5 mL/min	0.01 5 mL/min
カラム内径	L	0.5	1 mm	2.1 mm		3.0 mm	l	4.6 mm	

DAD 用のフローセルの選択

図 51 HPLC のフローセルの選択

フローセル光路長

Lambert-Beer の法則は、フローセルの光路長と吸光度が直線関係にあることを示しています。

Absorbance =
$$-\log T = \log \frac{I_0}{I} = \varepsilon \cdot C \cdot d$$

変数の意味は次のとおりです。

- ▼ はトランスミッションです。透過光線強度 I を入射光線強度 I₀ で割った 商として定義されたものです。
- e は吸光係数です。波長、溶媒、温度およびその他のパラメータが正確に定 義された条件下での、各々の物質の特性です。
- **C** は吸光試料の物質量(通常、g/L または mg/L 単位)。
- **d** は測定に使用するセルの光路長です。

適切なカラムとフローセルの組合せ

フローセルの光路長が長いほど、シグナルは強くなります。通常、光路長 が増加してもノイズは余り増加しないので、シグナル/ノイズ比が大きく なります。たとえば、ノイズが 10 % 未満増加するのに対し、シグナル強 度が 70 % 増加するのは、光路長が 6 mm から 10 mm に増加した場合で す。

光路長を増加させると、通常はセル容量が増加します。この例では、 5 μ L ~ 13 μ L の範囲です。通常、セル容量が増加するとピーク分散が大 きくなります。『142 ページ 図 52』に示すように、この例では、グラジェ ント分離の分解能に影響は及ぼしません。

目安として、フローセル容量は、半値幅におけるピーク量の約 1/3 である ことが理想です。ピーク量を計算するには、積分結果としてレポートされ たピーク幅に流速を掛け、その値を 3 で割ります。



図 52 シグナルの高さに及ぼすセル光路長の影響

従来、UV 検出器による LC 分析は、内部または外部標準と測定値の比較に 基づくものでした。検出器の測光精度をチェックするために、検出器フ ローセルの光路長に関するより正確な情報が必要となります。

正しいレスポンスは以下のとおりです。

予測レスポンス * 補正ファクタ

フローセルの詳細を 『143 ページ 表 24』と 『144 ページ 表 25』に示します。

適切なカラムとフローセルの組合せ

VWD フローセルの補正ファクタ

表 24 Agilent VWD フローセルの補正ファクタ

フローセルタイプ	セル容量	部品番号	光路長 (公称)	光路長 (実測値)	補正フ アクタ
標準フローセル	14 µL	標準フローセ ル、10 mm、 14 µL、40 bar (G1314-60086)	10 mm	10.15 ± 0.19 mm	10 /10.15
セミマイクロ フローセル	5 µL	セミマイクロ フローセル 6 mm、5 μL、 40 bar (G1314-60083)	6 mm	$6.10 \pm 0.19 \text{ mm}$	6 /6.10
マイクロ フローセル	2 μL	マイクロフ ローセル、 ID タグなし、 3 mm、2 µL、 120 bar (G1314-60087)	3 mm	2.80 ± 0.19 mm	3 /2.8
高圧フローセル	14 µL	高耐圧フロー セル 10 mm、 14 µL、400 bar (G1314-60082)	10 mm	10.00 ± 0.19 mm	6 /5.75

注記

機械加工誤差と比べると非常に小さいと思えるものですが、ガスケットの厚み とその圧縮率には誤差があることを意識する必要があります。

適切なカラムとフローセルの組合せ

DAD フローセルの補正ファクタ

表 25 フローセルの補正ファクタ

フローセル	光路長(実測値)	補正ファクタ
標準フローセル、10 mm、13 μL、120 bar (12 MPa) (G1315-60022)	9.80 \pm 0.07 mm	10/9.8
セミマイクロフローセル、6 mm、5 μL、120 bar (12 MPa) (G1315-60025)	5.80 \pm 0.07 mm	6/5.8
マイクロフローセル、3 mm、2 μL、120 bar (12 MPa) (G1315-60024)	3.00 +0.05 mm/-0.07 mm	3/3
セミナノフローセルキット、10 mm、500 nL、5 MPa (G1315-68724)	$10.00 \pm 0.02 \text{ mm}$	10/10
ナノフローセルキット、6 mm, 80 nL、5 MPa (G1315-68716)	$6.00 \pm 0.02 \text{ mm}$	6/6
標準フロー セルバイオー不活性、10 mm, 13 µL、 120 bar(12 MPa)MWD/DAD 用、キャピラリキットフ ローセル BIO(p/n G5615-68755)(G5615-60022)	9.80 \pm 0.07 mm	10/9.8

注記

機械加工誤差と比べると非常に小さいと思えるものですが、ガスケットの厚み とその圧縮率には誤差があることを意識する必要があります。
1220 Infinity LC

9

テスト機能とキャリブレーション

Agilent 1220 Infinity LC システム 147 据え付けチェック 147 モジュール情報 148 ステータス情報 148 溶媒送液システム 149 アイソクラティック ポンプのリークテストの説明 149 グラジエントポンプのリーク テストの説明 150 リーク テストの実行 152 リーク テスト結果の評価 153 高すぎる圧力のチェック 157 高すぎる圧力のチェックの評価 158 パージ ポンプ 159 オート サンプラ 160 メンテナンス ポジション 160 インジェクタの手順 161 アライメントティーチング 164 グリッパのベリフィケーション 167 カラムオーブン 168 オーブン テスト 168 オーブン キャリブレーション 169



Agilent Technologies

適切なカラムとフローセルの組合せ

可変波長検出器 (VWD) 170 セルテスト 170 暗雷流テスト 171 ホルミウムオキサイドテスト 172 強度テスト 173 フィルタ / グレーティング モーターテスト 176 検出器のキャリブレーション 177 クロマトグラムのテスト 178 スペクトルスキャン 179 ダイオードアレイ検出器 (DAD) 180 セルフテスト 180 フィルタ テスト 182 スリットテスト 184 暗電流テスト 185 強度テスト 187 ホルミウムオキサイドテスト 190 スペクトル平坦性テスト 193 ASTM ノイズテスト 194 セルテスト 195 内蔵テストクロマトグラムの使用 197 波長ベリフィケーションおよびリキャリブレーショ ン 199 Agilent ChemStation における診断情報 200 D/A コンバータ (DAC) テスト 202

この章では、Instrument Utilities ソフトウェアまたは Lab Advisor で利用できるテスト、キャリブレーションおよびツー ルについて説明します。

Agilent 1220 Infinity LC システム

Agilent 1220 Infinity LC システム

据え付けチェック

据え付けチェックでは、設定できるモジュールすべてをオンにし、システ ムを 1 mL/min で 5 分間パージし、最大 200 bar の圧力を加えて流路を テストし、オーブン(利用可能な場合)と検出器をオンにします。

注記 ポンプと検出器はこのチェックを必ず行う必要があります。オーブンとオート サンプラについては任意です。

以下の条件を満たしている場合、据え付けチェックは合格となります。

- すべてのモジュールがタイムアウト期間(120 s)内に正常にオンになる。
- ポンプが、5 min 後に 200 bar に達する。
- オーブンが実際より2K高い温度に達する。
- ・ 検出器ランプが点火し、検出器が レディ 状態に達する。

Agilent 1220 Infinity LC システム

モジュール情報

[モジュール情報]ツールは、モジュールからの診断情報を収集し、結果 をファイルに書き込みます。結果は次の3つのタブで表示できます。

一般

[一般] タブには、2 列の表形式で、モジュールのファームウェアとオプ ションに関する情報が表示されます。

テーブル

[テーブル] タブでは、そのモジュールに対応しているすべての診断テー ブルの内容が表示されます。[+] 記号をクリックするとテーブルが開き、 [-] 記号をクリックすると開いていたテーブルが閉じます。

シグナル

[シグナル] タブには、モジュールからの使用可能な診断シグナルのプ ロットが表示されます。使用可能なシグナルプロットは、モジュールごと に異なります。使用可能な場合は、1 つのシグナルにつき、短期的なプ ロットと長期的なプロットの両方が表示されます。

ステータス情報

[ステータス情報] ツールは、すべての Agilent 1220 Infinity LC モジュールの現在の状態を表示します。状態は継続的に更新されます。中断しない場合、60 分間、実行します。ツールを停止するには、[中断]をクリックします。

テスト機能とキャリブレーション 9 溶媒送液システム

溶媒送液システム

アイソクラティック ポンプのリークテストの説明

リークテストは、ポンプの耐圧性を確認するための内蔵トラブルシュー ティングテストです。テストには、あらかじめ定義したポンプシーケンス に従ってポンプを稼働したときの圧力プロファイルのモニタリングを行い ます。結果の圧力プロファイルから、ポンプコンポーネントの耐圧性と動 作に関する情報を得ることができます。

1段目:

初期化後、プランジャ 2 はストロークの最上位にあります。テストはプランジャ 1 から開始され、テスト時のストローク長は 100 μ L、流量は 153 μ L/min です。昇圧中のプランジャシーケンスは、1 -2 -1 -2 になります。

水平部 1:

プランジャ 2 は、約 1 分間、流量 2 μL/min でポンプに送液を続けま す。

2 段目:

流量は 153 µL/min に変更され、プランジャ 2 は残りのストローク分を 送液します。その後、プランジャ 1 が 2 段目の後半部分が完了するまで 送液を続けます。

水平部 2:

約 1 分間、流量が 2 µL/min に減らされます (プランジャ 1 が引き続き 送液)。 9 テスト機能とキャリブレーション 溶媒送液システム

3段目:

流量が 220 μL/min に増やされ、ストロークは 100 μL に変更されます。 プランジャ 1 がストロークを完了します。次に、流量が 510 μL/min に 変更されます。プランジャシーケンスは 2 -1 -2 -1 で、390 bar まで昇 圧されます。

水平部 3:

システム圧力が 390 bar に達した時点で、流量は 0 に変更し、圧力は 400 bar で安定させます。

最大圧力に達した後 1 分間経過しても、圧力低下が 2bar/min を超えてはいけません。

グラジエントポンプのリーク テストの説明

リークテストは、ポンプコンポーネントの耐圧性を確認するための内蔵ト ラブルシューティングテストです。テストは、ポンプで問題が発生したと 疑われる場合に実行します。テストでは、さまざまなプランジャが溶媒を 送液している間、非常に低流量で圧力増加をモニタリングします。あらか じめ設定された一連のポンプシーケンスを実行して、低流量における圧力 プロファイルを評価することにより、非常に小さなリークも検出できる。 テストでは、ポンプをブランクナットで詰まらせる必要があります。その 後、イソプロパノール(IPA)を使用してテストを実行し、同時に圧力プロ ファイルをモニタリングします。

注記 システムの加圧を始める前に、流路全体が IPA によって完全に洗浄されてい ることを必ず確認してください。流路中に、他の溶媒が残っていたり、微少な 気泡があると、テストは必ず失敗します。

溶媒送液システム

1段目

初期化後、プランジャ 2 はストロークの最上位にあります。テストはプ ランジャ 1 から開始され、テスト時のストローク長は 100 µL、流量は 153 µL/min です。昇圧中のプランジャシーケンスは、1 -2 -1 -2 になり ます。本来なら、この段階で圧力増加は直線的になるはずです。この段階 での圧力障害は、大きなリークまたはポンプコンポーネントの欠陥がある ことを示しています。

水平部1

プランジャ 2 は、約 1 分間、流量 2 μL/min でポンプに送液を続けま す。水平部での圧力は、一定の状態を保つか、わずかに増加となる必要が あります。圧力の低下は、2 μL/min を超えるリークを示しています。

2段目

流量は 153 μL/min に変更され、プランジャ 2 は残りのストローク分を 送液します。その後、プランジャ 1 が 2 段目の後半部分が完了するまで 送液を続けます。

水平部2

約 1 分間、流量が 2 µL/min に減らされます(プランジャ 1 が引き続き 送液)。水平部での圧力は、一定の状態を保つか、わずかに増加となる必要 があります。圧力の低下は、2 µL/min を超えるリークを示しています。

3段目

流量が 220 μ L/min に増やされ、ストロークは 100 μ L に変更されます。 プランジャ 1 がストロークを完了します。次に、流量が 510 μ L/min に 変更されます。プランジャシーケンスは 2 -1 -2 -1 で、390 bar まで昇 圧されます。

水平部3

システム圧力が 390 bar に達した時点で、流量は 0 に変更し、圧力は 400 bar で安定させます。

最大圧力に達した後 1 分間経過しても、圧力低下が 2 bar/min を超える ことはありません。 9 テスト機能とキャリブレーション 溶媒送液システム

リーク テストの実行

- **日時**: ポンプの問題が疑われる場合。
- **必要なツール: 説明** スパナ、1/4 インチ(HPLC ツールキットに付属)
- 必要な部品: 番号 部品番号 説明
 1 G1313-87305 リストリクションキャピラリ
 1 01080-83202 ブランクナット
 1 500 mL イソプロパノール
- **必要な準備**: ・ 溶媒キャビネットに、LC グレードのイソプロピルアルコールのボトルを セットして、その溶媒チューブをポンプのアクティブインレットバルブに 接続します。
- 注記 システムの加圧開始前に、テストする流路のすべての部品が IPA で完全にフ ラッシュされているか絶対に確認します。流路中に他の溶媒が残っていたり、 微少な気泡があると、テストは必ず失敗します。

Agilent Lab Advisor からのテスト実行

- 1 [テスト選択] メニューから [リークテスト] を選択します。
- 2 テストを開始して、指示に従います。
- 注記 テストが完了した際に、パージバルブをゆっくりと開けて、圧力を必ず開放し ます。
- 注記 『「リーク テスト結果の評価」153 ページ』で、リーク テスト結果の評価と解 釈方法を説明します。
- 注記 詳細は、Agilent Lab Advisor ツールを参照してください。

溶媒送液システム

リーク テスト結果の評価

ポンプヘッド内のコンポーネントに欠陥またはリークがあると、リークテ スト圧力プロットが変化します。代表的な障害モードについて、以下で説 明します。

注記 テスト中のエラーとテストの失敗との違いに注意してください。エラーとは、 テストの実行中に異常終了したことを意味しています。テストの失敗とは、テ スト結果が指定されたリミット値内ではなかったことを示しています。

注記

テストの失敗は、ブランクナット自体の損傷(締めすぎによる変形)に起因していることがよくあります。他の可能性のある原因を調べる前に、使用しているブランクナットが正常で、適度に締められていることを確認してください。

表 26 1 段目で圧力が増加しな	εl	,	١
-------------------	----	---	---

考えられる原因	修正措置
ポンプが稼働していない。	ログブックでエラーメッセージを確 認します。
パージバルブが開いている。	パージバルブを閉じて、テストを再 実行します。
継ぎ手の緩み、またはリーク。	すべての継ぎ手が締まっていること を確認するか、キャピラリを交換し ます。
溶媒ラインの接続が間違っている。	デガッサからの溶媒ラインが正しく 接続されていることを確認します。
パージバルブが汚れている。	パージバルブを開閉し、汚れを洗浄 します。洗浄後もリークする場合は、 バルブを交換します。
ポンプシールで大量(目で確認でき るほど)のリークがある。	ポンプシールを交換します。
アクティブインレットバルブ、アウ トレットバルブ、またはパージバル ブで大量(目で確認できるほど)の リークがある。	リークのあるコンポーネントがしっ かり取り付けられていることを確認 します。必要であればコンポーネン トを交換します。

- 9 テスト機能とキャリブレーション 溶媒送液システム
 - 表 27 Eカリミットに達していないが、水平部が水平またはポジティブ である

考えられる原因	修正措置
デガッサとポンプが十分に洗浄され ていない(ポンプヘッド内の空気)。	加圧下でデガッサとポンプをイソプ ロパノールを使用してパージします (リストリクションキャピラリを使 用)。
溶媒が適切でない。	イソプロパノールを取り付けます。 デガッサとポンプをパージします。

表	28	すべての水平部がネガー	ティ	ィブ	であ	る
~				-	$\sim \omega_{2}$	-

考えられる原因	修正措置
継ぎ手の緩み、またはリーク。	すべての継ぎ手が締まっていること を確認するか、キャピラリを交換し ます。
パージバルブの緩み。	パージバルブを締めます(14 mm ス パナ)。
パージバルブが汚れている。	パージバルブを開閉し、汚れを洗浄 します。洗浄後もリークする場合は、 バルブを交換します。
ポンプヘッドネジの緩み。	ポンプヘッドネジが締まっているこ とを確認します。
シールがリークしているか、プラン ジャに傷が付いている。	ポンプシールを交換します。プラン ジャに傷がないか確認します。傷が ある場合は交換します。
アウトレットバルブにリークがある。	アウトレットバルブを交換します。
ダンパーにリークがある。	ダンパーを交換します。

考えられる原因	修正措置
ポンプ内に空気があるか、新しい シールがまだ固定されていない。	加圧下でイソプロパノールを使用して ポンプを洗浄します(リストリクショ ンキャピラリを使用)。
アクティブインレットバルブの緩 み。	アクティブインレットバルブを締めま す (14 mm スパナ)。締めすぎないで ください。
ポンプヘッドネジの緩み。	ポンプヘッドネジが締まっていること を確認します。
アウトレットバルブの緩み。	アウトレットバルブ内のフィルタが正 しく取り付けられていることを確認し ます。アウトレットバルブを締めま す。
シールがリークしているか、プラ ンジャに傷が付いている。	ポンプシールを交換します。プラン ジャに傷がないか確認します。傷があ る場合は交換します。
アクティブインレットバルブの欠 陥。	アクティブインレットバルブを交換し ます。

表 29 1 番目の水平部がポジティブで、2 番目と 3 番目の水平部がネガ ティブである

1220 Infinity LC

9 テスト機能とキャリブレーション 溶媒送液システム

考えられる原因	修正措置
アウトレットバルブにリークがあ る。	アウトレットバルブをクリーニングし ます。アウトレットバルブ内のフィル タが正しく取り付けられていることを 確認します。アウトレットバルブを締 めます。
ポンプヘッドネジの緩み。	ポンプヘッドネジが締まっていること を確認します。
シールがリークしているか、プラ ンジャに傷が付いている。	ポンプシールを交換します。プラン ジャに傷がないか確認します。傷があ る場合は交換します。

表 30 1 番目の水平部がネガティブで、2 番目の水平部がポジティブで ある

表 31 3 段目がリミットに達しない

考えられる原因	修正措置
エラーが原因でポンプが停止した。	ログブックでエラーメッセージを確 認します。
ポンプシールで大量(目で確認でき るほど)のリークがある。	ポンプシールを交換します。
アクティブインレットバルブ、アウ トレットバルブ、またはパージバル ブで大量(目で確認できるほど)の リークがある。	リークのあるコンポーネントがしっ かり取り付けられていることを確認 します。必要であればコンポーネン トを交換します。

考えられる原因	修正措置
継ぎ手の緩み、またはリーク。	すべての継ぎ手が締まっていること を確認するか、キャピラリを交換し ます。
パージバルブの緩み。	パージバルブを締めます(14 mm ス パナ)。
パージバルブが汚れている。	パージバルブを開閉し、汚れを洗浄 します。洗浄後もリークする場合は、 バルブを交換します。
ポンプヘッドネジの緩み。	ポンプヘッドネジが締まっているこ とを確認します。
シールがリークしているか、プラン ジャに傷が付いている。	ポンプシールを交換します。プラン ジャに傷がないか確認します。傷が ある場合は交換します。
アウトレットバルブにリークがある。	アウトレットバルブを交換します。
ダンパーにリークがある。	ダンパーを交換します。

表 32 3 番目の水平部がネガティブである(圧力降下が 2 bar/min を超 えている)

高すぎる圧力のチェック

このテストでは、流路の詰まりをチェックし、詰まりの原因になっている モジュールの特定を試みます。詰まりがオートサンプラにある場合、テス トで詰まりがニードルで発生しているのかニードルシートで発生している のかを特定できます。

[高圧異常チェック]を実行するには、ポンプとオートサンプラが必要です。

高すぎる圧力のチェックの評価

開始条件

ポンプとオートサンプラをレディ状態にして、システムに動作圧力 200 bar を加えます。

ポンプリップルを測定し、リップルが定義されたリミット値(通常、動作 圧力の1%)内に入るまでテストの開始を遅らせます。

テスト第1部

テストの第1 部では、システムのどの部分で圧力の問題が発生している かの特定します。

システムで開始条件が揃った後、オートサンプラのバルブがメインパスか らバイパスに切り替えられ、圧力スライスがリミット値に対してテストさ れます。

リミット値を超える場合、圧力の問題はオートサンプラにあり、それ以外 の場合は、流路のどこかで問題が発生しています。

テスト第2部

テストの第2 部では、空のバイアルをニードルに向けて動かし、バルブ をバイパスからメインパスに切り替えて、圧力を大幅に低下させます。圧 力の低下をリミット値に対してチェックします。

オートサンプラが問題の原因と特定され、リミット値に達しない場合、問 題はニードル、ニードルループまたはメタリングドライブにあり、それ以 外の場合は、ニードルシートまたはニードルシートキャピラリにあります。

オートサンプラが問題の原因ではない場合、問題はポンプ(通常、フィル タまたはフリット)またはオートサンプラの後(ヒーターキャピラリまた はカラム)にあります。圧力下降のリミット値に到達しない場合、問題は ポンプはあり、それ以外の場合は、問題はオートサンプラの後で発生して います。

溶媒送液システム

パージ ポンプ

パージポンプの説明

[パージポンプ] ツールを使用すると、指定した時間に指定した流量でポンプから溶媒をパージできます。マルチチャンネルポンプ、および溶媒選択バルブ (SSV) 付きのポンプでは、パージするチャンネルを選択します。 各チャンネルは異なる条件でパージできます。

流量は、1 ~ 5 mL/min の範囲で 1 mL/min 刻みで選択できます。

時間は、1、2、3、5、7、10、15 分から選択できます。

G1361A 分取ポンプにはオートパージサイクルが用意されているため、ユー ザーが設定可能なオプションはありません。

ポンプのパージ

ポンプをパージするには

- 1 各チャンネルに適切なパージ溶媒を準備します。
- **2 [ツール選択]**画面で [パージポンプ]を選択します。
- **3** [パージ設定] ダイアログボックスで、
 - a 必要に応じて、パージするチャンネルを選択します。
 - **b** 選択したチャンネルごとに、[流量]およびパージする[時間]を選 択します。
 - c [OK] をクリックして、[パージ設定] ダイアログボックスを閉じま す。
- 4 パージバルブを開くようにリクエストが表示されたら、ポンプのパージ バルブを開き、[OK] をクリックして、メッセージボックスを閉じます。 パージ中、[一般] タブにパージされている現在のチャンネルと残って いるパージ時間が表示されます。[シグナル] タブには、パージサイク ル全体に対する時間対圧力のプロットが表示されます。
- 5 パージ時間が経過し、パージバルブを閉じるようにリクエストが表示されたら、ポンプのパージバルブを閉じ、[OK] をクリックして、メッセージボックスを閉じます。 これでポンプのパージプロセスが完了しました。

注記

9 テスト機能とキャリブレーション オート サンプラ

オート サンプラ

メンテナンス ポジション

ニードル交換

[**ニードル交換**]機能は、セーフティフラップを別の位置に移動し、ニードルを容易に交換およびアライメントできる位置に配置します。

[開始]	セーフティフラップをニードルから 離れた位置に移動し、ニードルを ニードルシートの約 15 mm 上に位置 付けます。
[アップ]	ニードルアームを徐々に上へ移動 します。
[ダウン]	ニードルアームを徐々に下へ移動し ます。ニードルをニードルアーム内 の正しい位置にアライメントするた めに、最も下の位置が使用されます。
[終了]	セーフティフラップをニードルを囲 む位置に再配置します。

ピストン交換

[ピストン交換]機能は、ピストンをホームポジションから離し、スプリングの張力を緩めます。この位置で、分析ヘッドアセンブリを取り外したり、メンテナンス後に簡単に再取り付けしたりすることができます。

[開始]	ピストンをホームポジションから離 し、スプリングの張力を緩めます。
[終了]	ピストンをホームポジションに戻し ます。

オート サンプラ

グリッパ交換

[**グリッパ交換**]機能は、グリッパをオートサンプラの前面に移動し、グリッパ開放メカニズムに簡単にアクセスできるようにします。

[開始]	グリッパをサンプルトレイ領域の前 面に移動します。
[終了]	グリッパをホームポジションに戻し ます。

アームポジション

[アームをホームへ移動]	トレイへのアクセスとトレイの交換 を容易にするため、グリッパアーム をホームポジションに移動します。
[アーム移動]	グリッパアームをサンプリングユ ニットの後ろのパーク位置に固定し ます。グリッパアームをパークする 前に、グリッパにバイアルがないこ とを確認してください。

インジェクタの手順

サンプリングシーケンスの各動作を手動コントロールで実行できます。こ れは、特定の障害モードを調べたり、修理が正常に完了したことを確認し たりするために、各サンプリングステップを緻密に観察する必要があるト ラブルシューティング時に役に立ちます。

各インジェクタステップコマンドは、実際には、一連の個々のコマンドから構成されています。個々のコマンドは、オートサンプラコンポーネントをあらかじめ定義した位置に移動し、特定のステップを実行できるようにします。

9 テスト機能とキャリブレーション オート サンプラ

表 33 インジェクタのステップコマンド

	.,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,	
ステップ	動作	注釈
[バルブバイパス]	注入バルブをバイパスの 位置に切り替えます。	
[プランジャホーム]	プランジャをホームポジ ションに移動します。	
[ニードルアップ]	ニードルアームを上部ポ ジションまで上げます。	バルブがバイパスポジションになっ ていない場合、バルブがバイパスに 切り換わります。
[バイアルをシートに移動]	選択したバイアルをシー トポジションまで移動し ます。	ニードルも上部のポジションに移動 します。
[ニードルをサンプルに挿入]	ニードルをサンプル内に 挿入します。	バイアルをバイアルをシートポジ ションに置き、ニードルを上部ポジ ションに移動します。
[吸引]	設定した注入量をメタリ ングデバイスが吸引しま す。	バイアルをシートポジションに置 き、ニードルを持ち上げ、ニードル をバイアルに挿入します。このコマ ンドは複数回実行できます(ただ し、最大吸引量が 100 µL を超える ことはできません)。メタリングデ バイスをリセットするには、[プラ ンジャホーム] を使用してくださ い。
[ニードルアップ]	ニードルをバイアルの外 に上げます。	バルブがバイパスポジションになっ ていない場合、バルブがバイパスに 切り換わります。
[バイアルをトレイに戻す]	選択したバイアルをトレ イポジションまで戻しま す。	ニードルも上部のポジションに移動 します。
[ニードルをシートに挿入]	ニードルアームをシート まで下ろします。	バイアルをトレイポジションまで戻 します。

162

テスト機能とキャリブレーション 9 オート サンプラ

表 33 インジェクタのステップコマンド

ステップ	動作	注釈
[バルブのメインパス]	注入バルブをメインパス ポジションに切り替えま す。	
[リセット]	インジェクタをリセット します。	

9 テスト機能とキャリブレーション オート サンプラ

アライメントティーチング

ALS アライメントティーチングの説明

アライメントティーチングツールは、修理のためにモジュールを分解した 後で発生する場合がある、オートサンプラグリッパ位置のわずかな偏差を 補正するために必要です。このツールを使用するには、100 バイアルトレ イをオートサンプラに挿入する必要があります。

アライメント手順では、リファレンスポイントとしてトレイの位置を 2 か 所使用します。トレイは長方形であり、2 つのポイントでアライメントす れば、トレイにある他のすべてのバイアル位置を修正できるためです。補 正計算が完了すると、X とシータの両方の値が小数点以下第 1 位に丸めら れます。アライメント手順の完了時に、修正されたグリッパ位置がモ ジュールのファームウェアに保存されます。

注記 オートサンプラの正しい操作を確保するには、アライメント手順を正しい順序 で完全に(どの部分もスキップせずに)実施する必要があります。

ボタン	説明	キーボードのショート カット
A	シータを増加してグリッ パを回転します。	[カーソル上]
◀	グリッパを水平に左に移 動します。	[カーソル左]
•	グリッパを水平に右に移 動します。	[カーソル右]。
▼	シータを減少してグリッ パを回転します。	[カーソル下]
[アームを上げる]	グリッパアームを上昇さ せます。	[Page Up]
[アームを下げる]	グリッパアームを下降さ せます。	[Page Down]

ALS アライメントコントロール

テスト機能とキャリブレーション 9 オート サンプラ

ボタン	説明	キーボードのショート カット
[グリッパを開く]	グリッパを開きます。	
[グリッパを閉じる]	グリッパを閉じます。	
[開始 <<]	手順の実行を開始しま す。起動時のみ表示され ます。	[Enter]
[続行 <<]	手順の次のステップに ジャンプします。アライ メント中のみ表示されま す。	[Enter]
[再実行]	ステップの実行を再開し ます。	

9 テスト機能とキャリブレーション オート サンプラ

ALS アライメントティーチングの実行

注記 オートサンプラの正しい操作を確保するには、アライメント手順を正しい順序 で完全に(どの部分もスキップせずに)実施する必要があります。

ALS をアライメントするには:

- 1 100 バイアルトレイをオートサンプラに挿入します。
- 2 位置 15 および 95 にキャップ付きバイアルを取り付けます。
- 3 [開始 <<] をクリックします。 グリッパアームが、バイアル 15 の上の位置に移動します。
- 4 補正値を工場出荷時のデフォルトにリセットするには、[はい]を、変 更しない場合は、[いいえ]を選択します。
- 5 [アームを下げる] を使用して、バイアル上部に接触しないようにグ リッパのフィンガをできる限り近付けます。
- 6 ▲ と ▼(回転)、および ◀ と ▶(左右の移動)を使用して、水平面でのグリッパの位置を調整します。
- **7 [グリッパを開く]**を使用して、グリッパのフィンガを開きます。
- 8 [アームを下げる]を使用して、バイアルのキャップがグリッパフィン ガのゴムと同じ高さになるまで、グリッパアームをさらに 5 mm 下げま す。
- 9 バイアルがグリッパフィンガの中心にあることを確認し、必要に応じて 位置を再調整します(ステップ 6)。
- 10 グリッパの位置が正しくなったら、[続行]をクリックします。
 - グリッパアームが、バイアル 95 の上の位置に移動します。
- 11 位置 95 でグリッパをアライメントするには、ステップ 6 ~ 9 を繰り 返します。
- 12 [続行]をクリックします。

計算が完了すると、X とシータの両方の値が小数点以下第 1 位に丸めら れます。補正値は、ここでサンプラの非揮発性メモリに保存され、サン プラが初期化されます。

テスト機能とキャリブレーション 9 オート サンプラ

グリッパのベリフィケーション

ALS グリッパベリフィケーションの説明

ベリフィケーション手順では、グリッパアライメントが正しいことを確認 するためのリファレンスポイントとして、いくつかのバイアル位置を使用 します。ベリフィケーションにより複数の位置がアライメント外である場 合は、アライメント手順を実行する必要があります。

グリッパポジションのベリフィケーション

バイアル位置 1 、10 、55 、81 、100 を位置ベリフィケーションに使用 できます。

- 1 確認対象の位置にあるバイアルトレイに空のキャップ付きバイアルを挿 入します。
- 2 [バイアル位置]メニューから1番目のバイアル位置を選択します。
- 3 [選択したポジションに移動]を選択します。
- 4 グリッパアームのフィンガがバイアルの上部中央にアライメントされている場合は、[バイアルを選択]ボタンを選択して、グリッパアームが正常にバイアルをトレイから外し上方に移動することを確認します。正確にアライメントされていない場合は、グリッパを再度アライメントする必要があります。
- 5 [バイアルを戻す] ボタンを選択して、グリッパがバイアルを正しく再 配置することを確認します。正確にアライメントされていない場合は、 グリッパを再度アライメントする必要があります。
- 6 次のバイアル位置に対して手順を繰り返します。

9 テスト機能とキャリブレーション カラムオーブン

カラムオーブン

オーブン テスト

オーブンテストの説明

オーブンテストは、2つのペルチェ要素の加熱性能を評価するために使用 されます。

開始温度から 10 K の範囲の加熱速度が測定されます。開始温度は、30 ℃ から 50 ℃ までの間で、次のように決定されます。

- 現在のオーブン温度が 30 ℃ より低い場合、オーブンは温度を 30 ℃ にしようとします。30 ℃ が開始温度になります。
- 現在のオーブン温度が 30 ℃ より高く 50 ℃ より低い場合、現在の オーブン温度が開始温度になります。
- 現在のオーブン温度が 50 ℃ を超えている場合、エラーメッセージが表示されます。テストを実施するには、オープンを 50 ℃ より低い温度に冷却する必要があります。

オーブンテストの評価

[オーブンテスト] 終了時に、左右のチャンネルの昇温のスロープが評価 されます。両方のスロープが 4 ℃/min 以上であればテストは合格です。

テスト機能とキャリブレーション 9 カラムオーブン

オーブン キャリブレーション

オーブンキャリブレーションの説明

オーブンキャリブレーションの手順では、オーブン温度をキャリブレーションされた外部測定デバイスと比較して測定できます。

通常、温度キャリブレーションは、機器の寿命中には不要ですが、地域の 規制基準を順守するために、2 ポイントキャリブレーションおよびベリ フィケーション手順を実行できます。

オーブンキャリブレーションの実行

注記

測定およびキャリブレーションの手順には、Heraeus Quat340 クォーツ表面温 度測定センサーなど、必要な分解能と精度を備えた測定デバイスを推奨してい ます。発注情報については、最寄の Agilent Technologies サポート窓口にお 問い合わせください。

- 1 キャリブレーションされた温度測定デバイスを取り付けます。
- ユーザーインターフェースで、[オーブンキャリブレーション]を選択 します。
- 3 オーブンが最初の設定値(40°C)に達するのを待ちます。
- 4 熱交換器の温度を測定し、値をフィールドに入力します。
- 5 オーブンが2番目の設定値(50°C)に達するのを待ちます。
- 6 熱交換器の温度を測定し、値をフィールドに入力します。
- 7 [OK] をクリックして、キャリブレーション値をオーブンに保存するか、
 [キャンセル] をクリックして、キャリブレーションプロセスを中断します。

9 テスト機能とキャリブレーション 可変波長検出器 (VWD)

可変波長検出器(VWD)

セル テスト

VWD セルテストの説明

[セルテスト]は、グレーティングがゼロオーダーポジションにあるとき に、サンプリングによって測定された重水素ランプの強度とリファレンス ダイオードによって測定された強度(フィルタ未処理および非対数)を比 較します。結果の強度比(サンプル対リファレンス)は、フローセルで吸 収された光線のアマウントの割合です。フローセルウィンドウの汚れおよ び汚染を確認するためにも、このテストを使用できます。テストが開始さ れると、ゲインは-1 に設定されます。溶媒の吸収による影響を排除する ために、フローセル内で水を使用してテストを実行する必要があります。

テストはマイクロフローセルを使用して実行しないでください。光強度が低減 注記 しテストが失敗することがあります。

VWD セルテスト結果の評価

強度比は、フローセルウィンドウの汚れ具合と、使用しているフローセル のタイプに依存します。比率が低ければ低いほど、光がフローセルで吸収 されている割合が大きいことを意味します。

表 34 考えられる原因は、フローセルの過度の吸光度です。

原因	修正作業
フローセルに吸光性溶媒または気泡 があります。	確実にフローセルが水で満たされ、 気泡がないようにします。
フローセルが汚れている。	フローセルウィンドウを交換する。

可変波長検出器(VWD)

暗電流テスト

WWD 暗電流テストの説明

暗電流テストでは、サンプル回路、およびリファレンス回路のリーク電流 を測定します。このテストは、非直線性または過剰なベースラインノイズ の原因となる、サンプルダイオード、リファレンスダイオード、あるいは A/D コンバータの回路の故障をチェックするために使用されます。テスト 中には、ランプのスイッチをオフにする(G1314A/B/C)か、光路中に シャッターを挿入します(G1314D/E/F)。次に、各ダイオードからのリーク 電流が測定されます。テスト結果は自動的に評価されます。

WWD 暗電流テストの評価

表 35

リミット値

サンプル回路	>7900	カウント
リファレンス回路	>7900	カウント

テストの失敗原因

表 36 サンプル回路のノイズがリミット値を超えている:

原因	修正措置
サンプルダイオードの欠陥	サンプルダイオードを交換します。
サンプル ADC ボードの欠陥。	サンプル ADC ボードを交換します。

表 37 リファレンス回路のノイズがリミット値を超えている:

原因	修正措置
リファレンスダイオードの欠陥。	リファレンスダイオードを交換しま す。
リファレンス ADC ボードの欠陥。	リファレンス ADC ボードを交換しま す。

ホルミウムオキサイドテスト

WWD ホルミウムオキサイドテストの説明

[ホルミウムオキサイドテスト]では、内蔵酸化ホルミウムフィルタの3 つの最大吸光度特性を使用して、波長精度を確認します(「波長キャリブ レーション」も参照してください)。テスト結果は自動的に評価され、酸化 ホルミウムフィルタのスペクトルが提供されます。溶媒の吸収による影響 を排除するために、フローセル内で水を使用してテストを実行する必要が あります。テストが完了すると、結果が自動的に表示されます。

ホルミウムオキサイドテスト レポート



VWD Holmium Test Results

							Specificat	ion	Measu	red	Result
Dev Dev	viation viation	from from	wavelength wavelength	1:	360.8	nm nm	-11 -11	nm nm	0.0	nm nm	Passed Passed
Dev	viation	from	wavelength	3:	536.4	nm	-11	nm	0.0	nm	Passed

可変波長検出器(VWD)

VWD ホルミウムオキサイドテストの評価

テストは機器ごとに評価され、測定された最大値が自動的に表示されます。 1 つ以上の最大値がリミット値から外れた場合は、テストは不合格です。

表 38 リミット値

最大吸光度	リミット値
360.8 nm	-1 \sim +1 nm
418.5 nm	$-1~\sim$ +1 nm
536.4 nm	-1 \sim +1 nm

強度テスト

VWD 強度テストの説明

[強度テスト] では、VWD 全波長範範囲(190 - 800 nm) にわたる UV ラ ンプの強度が測定されます。テスト結果は自動的に評価され、強度スペク トルが提供されます。テストでは、全波長範囲での最高強度、平均強度、 およ最低強度を評価します。このテストを実行して、ランプと光学系の性 能を判断します(『「VWD セルテストの説明」170ページ』も参照してくだ さい)。溶媒の吸収による影響を排除するために、フローセル内で水を使用 してテストを実行する必要があります。強度スペクトルの形は、主として ランプとグレーティングに依存します。したがって、強度スペクトルは機 器によってわずかに異なります。テストの完了時に、強度スペクトルと強 度値が表示されます。

注記 テストはマイクロフローセルを使用して実行しないでください。光強度が低減 しテストが失敗することがあります。

9 テスト機能とキャリブレーション 可変波長検出器 (VWD)

強度テスト レポート



VWD Intensity Test Results

specir		Measured	Result
Accumulated lamp on time Highest intensity > 3200 Average intensity > 1600 Lowest intensity > 64	000 cts 71 000 cts 9 000 cts 9	94.35 h 123680 cts 951488 cts 36384 cts	Passed Passed Passed

可変波長検出器(VWD)

WWD 強度テストの評価

表 39 範囲

ランプ強度	リミット値(カウント)
最大	>320000
平均	>160000
最小	>6400

表 40 テスト不合格の考えられる原因

原因	修正作業
ランプ オフ	ランプの電源を入れます。
ランプが古い。	ランプを交換します。
フローセルに吸光性溶媒または気泡 があります。	確実にフローセルが水で満たされ、 気泡がないようにします。
フローセルが汚れている。	『「VWD セルテストの説明」170 ペー ジ』を実行します。テストが不合格 の場合は、フローセルウィンドウを 交換します。

可変波長検出器(VWD)

フィルタ / グレーティング モーターテスト

VWD フィルタ / グレーティングテストの説明

フィルタモーターとグレーティングモーターの実際の位置は、リファレンス(センサー)位置からのステップ数として定義されます。[VWD フィルタ/グレーティングテスト]は、フィルタモーターおよびグレーティングモーターがリファレンス(センサー)位置に戻るために必要なモーターのステップ数をカウントします。リファレンス位置に到達するまでに必要なステップ数が予測されたステップ数と同じ場合は、テストは合格です。モーターが移動できなかった場合や、モーターのステップを見失った場合は、テストは失敗です。テスト結果は自動的に評価されます。

VWD フィルタ / グレーティングテストの結果

テスト不合格の考えられる原因

表 41 フィルタモータテスト

原因	修正作業
フィルタモータアセンブリの不良。	フィルタモータアセンブリを交換し ます。
VWM ボードの不良。	VWM ボードを交換します。

表 42 グレーティングモータテスト

原因	修正作業
フィルタモータアセンブリの不良。	フィルタモータアセンブリを交換し ます。
VWM ボードの不良。	VWM ボードを交換します。

テスト機能とキャリブレーション 9 可変波長検出器(VWD)

検出器のキャリブレーション

波長ベリフィケーション / キャリブレーション

検出器の [波長キャリブレーション]は、重水素のゼロオーダーポジショ ンと 656 nm (アルファエミッションライン) およびベータエミッションラ インで 486 nm のエミッションラインポジションを使用して行われます。 キャリブレーション作業には3つのステップを必要とします。まず、ゼロ オーダーポジションでグレーティングがキャリブレーションされます。ゼ ロオーダー最大値が検出されるステッピングモータのステップポジション が検出器に保存されます。次に、グレーティングが 656 nm での重水素エ ミッションラインに対してキャリブレーションされ、最大になるモーター ポジションが検出器に保存されます。最後に、グレーティングが 486 nm での重水素エミッションラインに対してキャリブレーションされ、最大に なるモーターポジションが検出器に保存されます。

波長ベリフィケーション / キャリブレーションには約 2.5 分 を要し、初期ド 注記 リフトにより測定が歪められるため、ランプ点灯後の最初の 10 分以内は無効 にされます。

ランプが ON になった際、重水素ランプの 656 nm エミッションラインが 自動的に検出されます。

検出器のキャリブレーション実施時期

検出器は工場でキャリブレーションされており、通常の使用条件ではリ キャリブレーションは必要ないはずです。しかし、以下の場合はリキャリ ブレーションした方が良いと思います。

- メンテナンス後(フローセルまたはランプ)
- 光学ユニット内のコンポーネントの修理後
- 光学ユニットまたは VWM ボードの交換後
- 年に1 度以上の定期的に(稼働時適格性確認/性能確認作業の前など)
- クロマトグラフの結果から、検出器のリキャリブレーションが必要なこ とを示した場合

9 テスト機能とキャリブレーション 可変波長検出器 (VWD)

クロマトグラムのテスト

VWM ボードに内蔵されているあらかじめ定義された内蔵テストクロマトグ ラムは、ダイオードからの通常のシグナルと同様に ADC を通じて処理さ れ、ADC とデータ処理システムのチェックに使用することができます。シ グナルは、アナログ出力と、GPIB 上で取得可能です。

注記

クロマトグラムの分析時間は、レスポンスタイム(ピーク幅)の設定値に依存 します。ストップタイムが設定されていない場合、クロマトグラムは連続して 繰り返します。

レスポンスタイム	ストップタイム
0.06 秒	0.8 分
0.12 秒	0.8 分
0.25 秒	0.8 分
0.50 秒	0.8 分
1.00 秒	1.6 分
2.00 秒	3.2 分(デフォルト設定)
4.00 秒	6.4 分
8.00 秒	12.8 分

テストクロマトグラムには 4 つのメインピークがあり、それぞれ次の吸光 度を持ちます。

ピーク	吸光度(概算)
1	38 mAU
2	100 mAU
3	290 mAU
4	20 mAU

可変波長検出器(VWD)

スペクトルスキャン

[スペクトルスキャン] ツールは、ダイオードアレイ検出器および可変波 長検出器 (DAD/MWD および VWD) で使用できます。指定した範囲の波長に わたりスペクトルをスキャンして、他のアプリケーション (Microsoft Excel など) で使用できる CSV (コンマ区切り値) 形式のファイルにデー タをエクスポートします。

スキャンパラメータ

- **UV ランプオン** UV ランプをオンにします。
- **ブランクスキャン** (VWD 専用) および [範囲の終了] フィールドで指定し、分離度は [ステップ] フィールドで指定し、分離度は [ステップ] フィールドで指定します。
- サンプルスキャン 指定した分離度で、指定した範囲の波長にわたりサン プルスペクトルをスキャンします。波長の範囲は、[範 囲の開始]および[範囲の終了]フィールドで指定し、 分離度は [ステップ]フィールドで指定します。
- **データのエクス** 他のアプリケーションで使用できるように、選択した ポート データを CSV 形式でエクスポートします。

9 テスト機能とキャリブレーション ダイオードアレイ検出器(DAD)

ダイオードアレイ検出器 (DAD)

この章では、検出器の内蔵テスト機能について説明します。

セルフ テスト

DAD セルフテスト(『181 ページ 図 53』を参照)により一連の個別テスト を実行し、結果を自動的に評価します。以下のテストが実行されます。

- ・ フィルタ テスト
- スリットテスト
- ・ 暗電流テスト
- 強度テスト
- 波長キャリブレーションテスト
- ホルミウムテスト
- スペクトル平坦性テスト
- ASTM ノイズテスト (オプション)

セルフ テストは、1 回または繰り返して実行できます。繰り返し実行する ように設定した場合、ユーザーが停止するまでテストは連続して順番に実 行します。間欠的に起こる問題のトラブルシューティングには、テストを 繰り返し実行することが役立ちます。

ASTM ノイズテストでは、検出器ベースラインノイズ(254 nm)を、 1 mL/min で水を送液しながら測定します。このテストには約 20 分要しま す。必要に応じてセルフ テストシーケンスに含めたり、除外することがで きます。

[セルフ テスト] ダイアログボックスでセルフ テストの設定が行われます。[シングルテスト] または [反復テスト] のいずれかを選択します。
 [ASTM ノイズテスト] チェックボックスをオンにすると、セルフ テストにノイズテストが含まれます。
ダイオードアレイ検出器(DAD)

Agilent G1315 Self Test Results

			More INFU
	Limits	Measured	Result
Filter test	0.0050.5 AU	0.27 AU	Passed
Slit test	0.71.3	1.08	Passed
Dark current test	012000 cts	79778026 cts	Passed
Min. intensity (190nm - 220nm)	> 2000 cts	18042 cts	Passed
Min. intensity (221nm - 350nm)	> 5000 cts	28451 cts	Passed
Min. intensity (351nm - 500nm)	> 2000 cts	16886 cts	Passed
Min. intensity (501nm - 950nm)	> 2000 cts	14683 cts	Passed
Max. intensity (190nm - 350nm)	< 450000 cts	89461 cts	Passed
Max. intensity (700nm - 950nm)	< 300000 cts	63609 cts	Passed
Max. intensity (D2 alpha line)	< 1200000 cts	169947 cts	Passed
Wavelength at 486.0nm	485.5486.5 nm	486.31 nm	Passed
Wavelength at 656.1nm	655.6656.6 nm	656.25 nm	Passed
Holmium test	-11 nm	0.30 nm	Passed
Spectral flatness	< 0.002 AU	0.0000 AU	Passed
ASTM noise (20 min. at 254nm)	<= 0.02 mAU	0.0096 mAU	Passed

図 53 セルフ テスト結果(レポート)

詳細は、以下のページの個々のテストを参照してください。

9 テスト機能とキャリブレーション ダイオードアレイ検出器(DAD)

フィルタ テスト

フィルタ テストでは、フィルタアセンブリの正しい動作を確認します。テ ストが開始されると、酸化ホルミウム フィルタが所定の位置に移動されま す。フィルタの移動中、吸光度シグナルがモニタリングされます。フィル タの端が光路を通過する時に、極大波長が測定されます。フィルタが所定 の位置に到達するとすぐに、(酸化ホルミウムの)極大波長が測定されま す。最後に、フィルタは光路の外に移動されます。移動中、フィルタの端 が光路を通過する時に、追加の極大波長が期待されます。(フィルタ移動 中に)フィルタアセンブリの端から生じる 2 つの最大値が測定され、ホル ミウムオキサイドの極大波長がリミット値内の場合、テストは合格です。

Instrument:	G1315C
Serial Number:	PP0000024
Operator:	Wolfgang
Date:	25.02.2005
Time:	14:55:42
File:	C:\CHEM32\2\DIAGNOSE\DAD_FILTER3.DGR

		Result Status
DAD Filter Test		
Date: 25.02.2005; Time: 14:55:42 Expected total time: approx 45s		
Expected total time. approx. 455.	i	
Test Procedure: 1. Measuring and evaluating filter test result		0.16 done

図 54 フィルタ テストの結果(レポート)

フィルタ テストの評価

フィルタ テスト不合格

ダイオードアレイ検出器(DAD)

Test Failed テスト不合格

考えられる原因

対策

- フィルターアセンブリ(レバーと フィルタアセンブリを取り付けます。 フィルター)が設置されていません。
- **2** フィルターモータの故障。 Agilent のサービス担当者に連絡し

てください。

Holmium Oxide Maximum out of Limits

ホルミウムオキサイドの最大値がリミット範囲外である。

考えられる原因

- ホルミウムオキサイドフィルタが 酸化ホルミウム フィルタを取り付け 取り付けられていない。
 ます。
- 2 フィルターの汚れ。 酸化ホルミウム フィルタを交換しま す。

ダイオードアレイ検出器(DAD)

スリットテスト

スリットテストでは、マイクロメカニカルスリットの正しい動作を確認し ます。テスト中、検出器がランプ強度の変動をモニタリングする間に、ス リットはすべてのスリット位置に移動されます。スリット位置が変更され る時に、強度低下(小さなスリットに移動)または強度上昇(大きなス リットに移動)が定義した範囲内になる必要があります。強度変化が期待 範囲外の場合、テストは失敗です。

スリットテスト評価

リミット値: 0.7 - 1.3

考えられる原因

- フローセルがまだ設置されてい フローセルを外します。
 る。
- 古いか Agilent 製以外のランプ 『「強度テスト」187ページ』を実行 します。ランプが古いか故障してい る場合、交換します。
- 3 スリットアセンブリの故障 スリットアセンブリを交換します。 (迷光)。
- **4** 検出器メインボードの故障。 検出器メインボードを交換します。
- 5 PDA/光学ユニットの故障。 光学ユニットを交換します。

ダイオードアレイ検出器 (DAD)

暗電流テスト

暗電流テストでは、各ダイオードからのリーク電流を測定します。このテ ストを使用して、特定波長で非線形を生じる恐れのあるダイオードの漏れ を確認します。テスト中、スリットアセンブリは暗電流位置に移動し、ダ イオードアレイに達する光を遮断します。次に、各ダイオードからのリー ク電流が測定され、グラフに表示されます(『185 ページ 図 55』を参照)。 各ダイオードのリーク電流(カウントで表示)は、プロットに示されたリ ミット値(赤色の帯)内に入るはずです(『185 ページ 図 55』を参照)。

暗電流テスト評価

リミット値: 0... 12,000 カウント

Instrument:	G1315C	
Serial Number:	PP0000024	
Operator:	Wolfgang	
Date:	25.02.2005	
Time:	14:50:12	
File:	C:\CHEM32\2\DIAGNOSE\DAD	DARK2.DGR

Dark Current Plot



Dark Current Test Results

	Specification	Measured	Result
Dark current maximum value	<= 12000 cts	8056 cts	Passed
Dark current minimum value	> 0 cts	7994 cts	Passed

図 55 暗電流テストの結果(レポート)

9 テスト機能とキャリブレーション ダイナードアレイ検出器 (DAD)

ダイオードアレイ検出器 (DAD)

Test Failed テスト不合格

考えられる原因

- スリットアセンブリの故障 (迷光)。
- 2 検出器メインボードの故障。
- **3** PDA/ 光学ユニットの故障。

対策

『「スリットテスト」184 ページ』 (『「セルフ テスト」180 ページ』の 一部)を実行します。

- Agilent のサービス担当者に連絡し てください。
- Agilent のサービス担当者に連絡し てください。

ダイオードアレイ検出器(DAD)

強度テスト

注記

このテストは標準フロー セル (10 mm および 6 mm の光路長) 専用です。ナノ フローセル (80 nL および 500 nL) は、低容量のためこのテストを行うこと ができません。

強度テストでは、重水素ランプとタングステンランプの強度を全波長範囲 (190 - 950 nm) にわたって測定します。4 つのスペクトル範囲を使用し て、強度スペクトルを評価します。このテストを実行して、ランプと光学 系の性能を判断します(『「セル テスト」195 ページ』も参照)。テストを 開始すると、1-nm スリットが光路に自動的に移動し、ゲインがゼロに設定 されます。吸収を持つ溶媒の影響を除去するために、テストはフローセル を水で満たして行います。強度スペクトルの形は、主としてランプ、グ レーティング、およびダイオードアレイの特性に依存します。そのため、 機器間で強度スペクトルが微妙に異なります。『188 ページ 図 56』には、 代表的な強度テストのスペクトルを示します。

強度テストの評価

Agilent Lab Advisor、ChemStation、インスタント パイロットは、4 つの スペクトル範囲について自動的に評価を行い、各範囲のリミット値、測定 された強度カウント、そして各スペクトル範囲に対して合格または不合格 を表示します(『188ページ図 56』を参照)。 9 テスト機能とキャリブレーション ダイオードアレイ検出器 (DAD)

Instrument:	G1315C
Serial Number:	DE60755000
Operator:	
Date:	3/17/2009
Time:	1:24:38 PM

Intensity Plot



Intensity Test Results

	Specification	Measured	Result
Accumulated UV lamp on time		68.08 h	
Lowest intensity in range 190nm - 220nm:	> 2000 cts	18069 cts	Passed
Lowest intensity in range 221nm - 350nm:	> 5000 cts	28464 cts	Passed
Lowest intensity in range 351nm - 500nm:	> 2000 cts	16889 cts	Passed
Lowest intensity in range 501nm - 950nm:	> 2000 cts	14665 cts	Passed
Highest intensity in range 190nm - 350nm	: < 450000 cts	89478 cts	Passed
Highest intensity in range 700mm - 950mm	: < 300000 cts	63598 cts	Passed
Highest intensity for the D2 alpha line:	< 1200000 cts	169933 cts	Passed

図 56 強度テストの結果 (レポート)

1 つ以上の範囲でカウントが低い場合は、フローセルを使用した場合とフ ローセルを取り除いた場合の値を比較するテストを開始します。

セル窓 / レンズ (Vis- ランプとフローセルの間に 3 個あります) が汚れ ていると、光のスループットは減少します。

ダイオードアレイ検出器(DAD)

もし検出器が、501 nm ~ 950 nm の範囲で失敗した場合は、以下を確認し てください。

- Vis-ランプはオンになっていますか? オンになっていなければ、オン にします。
- Vis-ランプのガラスが汚れているか、壊れていますか? 汚れたり壊れている場合には、Vis-ランプを交換します。
- UV- ランプは、Vis- ランプに面して反射コーティングを示しますか? 該 当する場合は、UV- ランプを交換します。
- 例 (フローセル無しで測定):

Vis- ランプ オフ、または不具合:110 カウントVis- ランプ オン、OK:13613 カウント

Test Failed テスト不合格

考えられる原因

- フローセルの中に吸光性溶媒また 確実にフローセルが水で満たされ、 は気泡が入っている。
 気泡がないようにします。
- 2 フローセルが汚れている。 セル テストを実行します(『「セル テスト」195ページ』を参照)。テストが不合格の場合は、フローセル窓 を交換します。
- 3 光学部品(アクロマート、ウィン 光学系のコンポーネントをアルコードウ)が汚れている。
 ルと糸くずが出ない布できれいにするか、または部品を交換します。
- **4** 古いか Agilent 製以外のランプ ランプを交換します。

ダイオードアレイ検出器(DAD)

注記 もしランプが、1 つの範囲で失敗した場合、アプリケーションをその特定の範 囲で実行しないのであれば、ランプを交換する必要はありません。

フローセルを取り外してテストをやり直します。もしカウントが劇的に上昇した場合(2 倍以上であれば、フローセルのコンポーネントは汚れており、メン テナンス / サービスが必要になることがあります。

ランプ交換の間隔が短くなっている場合、Agilent サービス担当が光学ユニットの光路にある汚れたコンポーネント(カップリング レンズ、ソース レンズ、セルサポート アセンブリ、フローセル窓)をチェックします。

ホルミウムオキサイドテスト

ホルミウムオキサイドテストでは、内蔵の酸化ホルミウム フィルタの 3 つの最大吸光度特性を使用して、波長真度を確認します(『「波長ベリ フィケーションおよびリキャリブレーション」199ページ』も参照)。テス トを開始すると、1 nm スリットが光路に自動的に移動します。吸収を持つ 溶媒の影響を除去するために、テストはフローセルを水で満たして行いま す。

『「ホルミウムフィルタの波長の証明書」417ページ』も参照してください。

ホルミウムオキサイドテストの評価

リミット値:

361.0 nm360.0 \sim 362.0 nm (± 1
nm)453.7 nm452.7 \sim 454.7 nm (± 1
nm)536.7 nm535.7 \sim 537.7 nm (± 1
nm)

注記

ダイオードアレイ検出器 (DAD)

テストは機器ごとに評価され、測定された最大値が自動的に表示されます。 1 つ以上の最大値がリミット値から外れた場合は、テストは不合格です (『191 ページ 図 57』を参照)。

Instrument:	G1315C
Serial Number:	PP0000024
Operator:	Wolfgang
Date:	25.02.2005
Time:	14:30:08
File:	C:\CHEM32\2\DIAGNOSE\DAD_FILTER2.DGR



Holmium Filter Test Results

		Specification	Measured	Result
Wavelength 1: 361.0	nm	360362 nm	360.9 nm	Passed
Wavelength 2: 453.7	nm	452.7454.7 nm	453.4 ni	m Passed
Wavelength 3: 536.7	nm	535.7537.7 nm	536.8 ni	m Passed

図 57 ホルミウムオキサイドテスト結果(レポート)

ダイオードアレイ検出器(DAD)

- Test Failed
- テスト不合格

考えられる原因

は気泡が入っている。

3 フローセルが汚れている。

- 1 フローセルの中に吸光性溶媒また フローセルが水で満たされているか を確認します。
- 2 キャリブレーションが正しくな リキャリブレーションを行い(『「波 長ベリフィケーションおよびリキャ い。
 - リブレーション」199ページ』を参 照)、テストを繰り返します。
 - セル テストを実行します (『「セル テスト」195ページ』を参照)。テス トが不合格の場合は、フローセル窓 を交換します。
- 4 光学部品(アクロマート、ウィン 光学系のコンポーネントをアルコー ルと糸くずが出ない布できれいにす ドウ)が汚れている。 るか、または部品を交換します (『「強度テスト」187ページ』を参 照)。
- 5 古いか Agilent 製以外のランプ UV ランプを交換します。

ダイオードアレイ検出器(DAD)

スペクトル平坦性テスト

スペクトル平坦性テストでは、スペクトルの最大ノイズ(単位 mAU)を測 定します。このテストはフローセルを取り外した状態で実行し、吸収を持 つ溶媒またはフローセルの汚れによる影響を排除します。まず、検出器の 調整を行います。次に、スペクトル(フローセルなし)を取得します。

理論的にはスペクトルは平坦ですが、実際にはノイズがスペクトル上に重 ね合わされます。ノイズの大きさは、光学系や電子機器の性能の指標です。

注記 このテストは検出器自己診断のみの一部です(『「セルフ テスト」180 ページ』 を参照)。

スペクトル平坦性評価

リミット値は、< 0.002 AU です。

考えられる原因

2 DAM ボードの欠陥。

- 1 古いか Agilent 製以外のランプ ・ 強度テストを実行します。
 - ランプが古いか故障している場合、交換します。
 - DAM ボードを交換します。
- 3 PDA/光学ユニットの故障。 光学ユニットを交換します。

9 テスト機能とキャリブレーション ダイオードアレイ検出器(DAD)

ASTM ノイズテスト

ASTM ノイズテストでは、20 分間にわたり検出器ノイズを測定します。フ ローセルを取り外した状態でテストを行うため、テスト結果は溶媒やポン プの影響を受けません。テストが完了すると、ノイズ結果が自動的に表示 されます。

注記

このテストは検出器自己診断のみの一部です(『「セルフ テスト」180ページ』 を参照)。

ASTM ノイズテストの評価

リミット値は ± 0.02 AU です。

考えられる原因

- ランプのウォームアップ時間が不 1時間以上、ウォームアップするた 十分。
 めにランプを放置します。
- **2** 古いか Agilent 製以外のランプ ランプを交換します。

ダイオードアレイ検出器(DAD)

セル テスト

セルテストでは、全波長範囲(190 ~ 950 nm)にわたる重水素ランプおよ びタングステンランプの強度を、フローセルが取り付けて 1 回、フローセ ルを取り外して 1 回測定します。結果の強度比は、フローセルで吸収され た光線のアマウントの割合です。フローセルウィンドウの汚染およびコン タミネーションを確認するためにも、このテストを使用できます。テスト を開始すると、1 nm スリットが光路に自動的に移動し、ゲインがゼロに設 定されます。吸光性溶媒の影響を除去するために、テストはフローセルを 水で満たして行います。

注記

新しい検出器 / フローセルには、最初にこのテストを行う必要があります。値は、今後の参照 / 比較のために保存する必要があります。

セル テストの評価

Agilent ChemStation により、強度比が自動的に計算されます。強度比 (新しい標準フロー セルで通常 $0.5 \sim 0.7$ 、新しいマイクロフローセルと 高圧フローセルで $0.1 \sim 0.3$)は、フローセルウィンドウの汚れ具合、お よび使用しているフローセルのタイプに依存します。

Instrument:	G1315C
Serial Number:	PP00000024
Operator:	Wolfgang
Date:	25.02.2005
Time:	14:54:22
File:	C:\CHEM32\2\DIAGNOSE\DAD_CELL2.DGR

	1	Result	Status
Detector Cell Test Date: 25.02.2005; Time: 14:54:22 Expected total time: approx. 45 s.			
Test Procedure: 1. If cell not in place, install it 2. Measuring intensity with cell 3. Remove cell 4. Measuring intensity without cell 5. Calculating intensity ratio		42053 42034 1.0	 done done done done done

図 58 フローセルを挿入しないセルテストの結果(レポート)

注記

標準フロー セルに対してだけ、このテストを使用できます。ナノフローセル では、その設計により非常に低い値になります。

ダイオードアレイ検出器(DAD)

テスト不合格(比の値が低すぎる)

考えられる原因

- フローセルに吸光性溶媒または気 確実にフローセルが水で満たされ、 泡があります。
 流行がないようにします。
- 2 フローセルが汚れている。 フローセルウィンドウを交換する。

ダイオードアレイ検出器(DAD)

内蔵テストクロマトグラムの使用

この機能は、Agilent ChemStation、Lab Advisor、インスタント パイロッ トから使用できます。

内蔵テストクロマトグラムを使用して、検出器からデータ システムやデー タ解析、またはアナログ出力を介したインテグレータやデータ システムへ のシグナルパスをチェックできます。クロマトグラムは、終了時間または マニュアル操作のいずれかにより終了を実行するまで連続的に繰り返しま す。

注記

ピーク高さは常に同じですが、面積とリテンションタイムは設定のピーク幅に より異なります。下の例を参照してください。

Agilent Lab Advisor を使用した手順

この手順は、すべての Agilent 1200 Infinity 検出器 (DAD、MWD、VWD、 FLD、RID) で行えます。例の図は、RID 検出器からのものです。

- 1 Assure デフォルト LC メソッドが、コントロール ソフトウェアにロー ドされていることを確認します。
- Agilent Lab Advisor ソフトウェア (B. 01. 03 SP4 以降)を開始し、検 出器のツール選択を開きます。
- **3** テストクロマトグラムの画面を開きます。

Tools: Test Chromatogram			
Current Status	Disabled	Switch Test Chromatogram on	
		Switch Test Chromatogram off	

- **4 テストクロマトグラム**をオンにします。
- 5 検出器のモジュール サービスセンター に変更し、検出器のシグナルを シグナル プロットウィンドウに追加します。

ダイオードアレイ検出器 (DAD)

6 テストクロマトグラムを開始するために、コマンド ラインに次を入力 します:STRT



7 テストクロマトグラムを終了するために、コマンド ラインに次を入力 します: STOP

注記

テストクロマトグラムは、分析終了時に自動的に電源が切れます。

ダイオードアレイ検出器(DAD)

波長ベリフィケーションおよびリキャリブレーション

検出器は、重水素ランプのアルファエミッション ライン(656.1 nm)と ベータエミッション ライン(486 nm)を使用して波長キャリブレーション を行います。シャープなエミッションラインによって、酸化ホルミウムを 使用する場合より正確なキャリブレーションが可能です。ベリフィケー ションを開始すると、1 nm スリットが光路に自動的に移動し、ゲインはゼ ロに設定されます。吸収を持つ溶媒の影響を除去するために、テストはフ ローセル内をバブルフリーの脱気した HPLC グレードの水で満たして行い ます。

偏差が確認および表示される場合、[調整]を押してリキャリブレーショ ンできます。偏差は [キャリブレーション履歴](検出器の診断バッファ) に記録されます。

	Devia D2-Alpha 0.01nm	Calibratio ation D2-Beta 0.04nm	on history Time	Date
			10.10.01	16.03.2005
D2-Alpha -1.2nm 656.1nm +1.2nm D2-Beta -1.2nm 486.0nm +1.2nm -1.2nm 486.0nm +1.2nm		Reset	history	
D2-Alpha line deviation is 0.00nm D2-Beta line deviation is 0.00nm Adjust	OK	Cancel	<u>H</u> elp	

図 60 波長ベリフィケーションおよびリキャリブレーション

次の作業後に、波長キャリブレーションを行う必要があります

- フローセルのメンテナンス後、
- ランプの交換、または
- プロセッサボードまたは光学ユニットの交換のような大きな修理の後、 『「モジュールのファームウェアの交換」361ページ』も参照してください。

キャリブレーションの後、ホルミウムオキサイドテスト(『191ページ図 57』を参照)により、3つの追加波長における波長真度の確認を行います。

Agilent ChemStation における診断情報

Agilent ChemStation は、検出器のさまざまなアセンブリに関する特定情報を提供します。

- 1 [診断]を開き、詳細表示に変更します。
- 2 対象アセンブリをクリックし、「変数表示の更新」を選択します。『201 ページ 図』に例を示します。

『200ページ表43』に詳細情報を示すアセンブリを一覧表示します。

表 43 診断 - 詳細情報

アセンブリ	詳細		
一般	製品番号、ファームウェアリビジョン、製造日 (機器、または、交換した場合はメインボードの)、 シリアル番号、積算点灯時間、分光計シリアル番 号。		
Vis ランプ	積算点灯時間、点灯時間実測値、ランプ点灯、積算 点灯時間スイッチ、電圧 / 電流のオン、ランプ電 流、ランプ電圧 / 電流		
UV ランプ	積算点灯時間、点灯時間実測値、ランプ点火、積算 点灯時間リセット、点火電圧 / 電流、ランプ電圧、 ランプ電圧 / 電流		
	ID タグ付きランプのみで使用可能。 製品番号、シリアル番号、製造日、最終強度テスト 日		
ホルミウムフィルター	フィルターの動き、フィルターの動きリセットログ		
フローセル	ID タグ付きフローセルのみで使用可能。 製品番号、製造日、光路長、最高圧力、シリアル番 号、容量、最終セルテスト日		
マイクロスリット	スリットの動き、フィルターの動きリセットログ		

ダイオードアレイ検出器 (DAD)



ダイオードアレイ検出器(DAD)

D/A コンバータ (DAC) テスト

検出器では、インテグレータ、チャートレコーダ、データ システムで使用 するために、クロマトグラフ シグナルのアナログ出力ができます。アナロ グ信号は、デジタルアナログコンバータ (DAC) によって、デジタル形式か ら変換されます。

DAC テストは、デジタルテスト シグナルを DAC に適用することによって、 デジタルアナログコンバータの正しい操作の確認に使用されます。

DAC は、約 50 mV のアナログ信号(もしアナログ出力のゼロオフセットが 初期値 5 % に設定されている場合)を出力します。これはインテグレータ にプロットできます。振幅 10 μ V および周波数約 1 cycle/24 seconds の 連続方形波がシグナルに適用されます。

矩形波の振幅とピーク - ピークノイズは、DAC テストの評価に使用されます。

- **日時**: アナログ検出器シグナルにノイズが多いか、不明な場合。
- **必要な準備:** ランプを 10 分以上オンにしておく必要があります。インテグレータ、 チャートレコーダ、またはデータシステムを検出器のアナログ出力に接続し ます。

Agilent LabAdvisor によるホルミウムオキサイドテストの実行

1 D/A コンバータ (DAC) テストを実行します (詳細については、ユー ザーインタフェースのオンラインヘルプを参照)。

Test N	ame	D/A Converter Test		Description	The test switches a test signal to the analog output, that can
Module	•	G4212A:PR00100015			be measured using an integrator or strip-chart recorder.
Status		Passed			
Start T	ime	7/9/2009 3:06:30 PM			
Stop T	ime	7/9/2009 3:06:53 PM			
Test Pro	cedure —				
💅 1.	Check P	rerequisites			
V 2.	Switch o	n Analog Output			
🖌 3.	Switch o	ff Analog Output			
図 62		D/A コンバータ	(DAC)	テスト - 糸	吉果

1220 Infinity LC

ダイオードアレイ検出器 (DAD)



インスタント パイロットによるフィルタモータテストの実行

コマンドラインにより、テストを開始できます。

- 1 テストを開始するには TEST: DAC 1 Reply: RA 00000 TEST:DAC 1
- 2 テストを停止するには TEST:DAC 0 Reply:RA 00000 TEST:DAC 0

Test Failed

テスト不合格

D/A コンバータ (DAC) テスト評価

このステップでのノイズは3 µV 未満である必要があります。

考えられる原因

対策

- 検出器と外部デバイス間のケーブ ケーブルをチェックするか取り替え ル不具合または接地に問題があ ます。 る。
- 検出器メインボードの故障。
 Agilent のサービス担当者に連絡してください。

1220 Infinity LC

9 テスト機能とキャリブレーション ダイオードアレイ検出器 (DAD)



エラー情報

10



エラー メッセージ内容 208 一般エラー メッセージ 209 Compensation Sensor Open 209 Compensation Sensor Short 210 Fan Failed 211 Leak 212 Leak Sensor Open 213 Leak Sensor Short 214 Remote Timeout 215 Shutdown 216 Lost CAN Partner 217 218 Timeout ポンプのエラー メッセージ 219 Encoder Missing 219 Index Adjustment 220 Index Limit 221 Index Missing 222 Initialization Failed 223 Missing Pressure Reading 223 Motor-Drive Power 224 225 Pressure Above Upper Limit Pressure Below Lower Limit 226 Pressure Signal Missing 227 Pump Configuration 227 Pump Head Missing 228 Restart Without Cover 228 Servo Restart Failed 229 Stroke Length 230



205

10 エラー情報 ダイオードアレイ検出器 (DAD)

> Temperature Limit Exceeded 231 Temperature Out of Range 232 Valve Failed (MCGV) 233 MCGV Fuse 234 Wait Timeout 235 Solvent Zero Counter 236 オート サンプラのエラー メッセージ 237 Arm Movement Failed 237 Initialization Failed 239 Initialization with Vial 240 Invalid Vial Position 241 Metering Home Failed 242 Missing Vial 243 Missing Wash Vial 244 Motor Temperature 245 Needle Down Failed 246 Needle Up Failed 247 Safety Flap Missing 248 Valve to Bypass Failed 249 Valve to Mainpass Failed 250 Vial in Gripper 251 一般検出器エラー メッセージ 252 252 Heater at fan assembly failed Heater Power At Limit 253 Illegal temperature value from sensor at fan assembly 253 Illegal Temperature Value from Sensor at Air Inlet 254 UV Lamp Current 255 UV Lamp Voltage 256

エラー情報 10 ダイオードアレイ検出器 (DAD)

VWD 検出器エラーメッセージ 257 ADC Hardware Error 257 Wavelength calibration setting failed 258 Cutoff filter doesn't decrease the light intensity at 250 nm 259 Filter Missing 260 Grating or Filter Motor Errors 261 Grating Missing 262 No heater current 263 Wavelength holmium check failed 264 Ignition Failed 265 Wavelength test failed 266 DAD 検出器エラー メッセージ 267 DAD 検出器エラー メッセージ 267 Diode Current Leakage 268 UV Ignition Failed 269 UV Heater Current 270 Calibration Values Invalid 271DAD 検出器エラー メッセージ 267 Wavelength Recalibration Lost 273 DSP Not Running 274 DAD 検出器エラー メッセージ 267 機器ログブック 275

この章では、表示されるエラーメッセージに関する情報を提供し、 考えられる原因とそれらを解決するための提案を示します。

10 エラー情報 エラー メッセージ内容

エラー メッセージ内容

分析を続けるために何らかの処置(修理、消耗品の交換など)を必要とす る障害が、電子部品、機械部品、および流路に発生した場合、ユーザーイ ンターフェイスにエラーメッセージが表示されます。このような障害が発 生した場合、モジュール前面の赤色ステータスインジケーターが点灯し、 モジュールログブックにエントリが書き込まれます。

エラー情報 10 一般エラー メッセージ

一般エラー メッセージ

一般エラーメッセージは、すべての Agilent シリーズ HPLC モジュールで 汎用的に使用されます。その他のモジュールでも同様に表示されることが あります。

Compensation Sensor Open

Error ID: 0081

補正センサーオープン

モジュールのメインボード上の周囲温度補正センサー (NTC)が故障しました (断線)。

メインボード上の温度補正センサー (NTC)の抵抗は、周囲温度によって 変化します。リーク回路は、この抵抗の変化を使用して、周囲温度の変化 を補正します。補正センサーの抵抗が上限値を超えた場合は、このエラー メッセージが生成されます。

考えられる原因

対策

1 メインボードの故障。 Agilent のサービス担当者に連絡し てください。



Compensation Sensor Short

Error ID: 0080

補正センサーショート

モジュールのメインボード上の周囲温度補正センサー(NTC)が故障しました (短絡)。

メインボード上の温度補正センサー (NTC)の抵抗は、周囲温度によって 変化します。リーク回路は、この抵抗の変化を使用して、周囲温度の変化 を補正します。センサーの抵抗が下限値を下回ると、このエラーメッセー ジが生成されます。

考えられる原因

対策

1 メインボードの故障。

Agilent のサービス担当者に連絡し てください。

エラー情報 10 一般エラー メッセージ

Fan Failed

Error ID: 0068

ファン動作不良

モジュールの冷却ファンが故障しました。

メインボードは、ファンシャフト上のホールセンサーを使用して、ファン の回転速度をモニタリングします。ファンの回転速度が一定期間、特定の リミット値以下に低下すると、エラーメッセージが生成されます。

モジュールによっては、アセンブリ(検出器内のランプなど)の電源がオ フとなることで、内部のモジュールが過熱するのを防ぎます。

考えられる原因

1	ファンケーブルの断線。	Agilent のサービス担当者に連絡し てください。
2	ファンの故障。	Agilent のサービス担当者に連絡し てください。
3	メインボードの故障。	Agilent のサービス担当者に連絡し てください。

Leak

Error ID: 0064

リーク

モジュールでリークが検出されました。

リークアルゴリズムが、2 つの温度センサー(リークセンサーとボード搭載の温度補正センサー)からのシグナルを使用して、リークが発生しているかどうか判断します。リークが発生すると、リークセンサーが溶媒によって冷却されます。これによるリークセンサーの抵抗の変化が、メインボード上のリークセンサー回路によって検知されます。

考えられる原因

1	フィッティングの緩み。	すべてのフィッティングがしっかり 締まっていることを確認します。
2	キャピラリの破損。	破損したキャピラリを交換します。

Leak Sensor Open

Error ID: 0083

リークセンサーオープン

モジュール内のリークセンサーが故障しました(オープン:断線)。

リークセンサーを流れる電流は、温度によって変化します。リークセン サーが溶媒によって冷却され、リークセンサー電流が規定のリミット値内 で変化したとき、リークが検出されます。リークセンサー電流が下限値よ り下がった場合は、このエラーメッセージが生成されます。

考えられる原因

1	リークセンサーがメインボードに 接続されていない。	Agilent のサービス担当者に連絡し てください。
2	リークセンサーの故障。	Agilent のサービス担当者に連絡し てください。
3	リークセンサーが正しく配線され ず、金属部品にはさまれている。	Agilent のサービス担当者に連絡し てください。



Leak Sensor Short

Error ID: 0082

リークセンサーショート

モジュールのリークセンサーが故障しました(短絡)。

リークセンサーを流れる電流は、温度によって変化します。リークセン サーが溶媒によって冷却され、リークセンサー電流が規定のリミット値内 で変化したとき、リークが検出されます。リークセンサー電流が上限値を 超えた場合は、このエラーメッセージが生成されます。

考えられる原因 対策 1 リークセンサーの故障。 Agilent のサービス担当者に連絡してください。 2 リークセンサーが正しく配線されず、金属部品にはさまれている。 Agilent のサービス担当者に連絡してください。

Remote Timeout

Error ID: 0070

リモートタイムアウト

リモート入力上にノットレディ状態が残っています。分析を開始すると、 通常は分析の開始から1分以内にすべてのノットレディ状態(検出器バラ ンス時など)がラン状態に切り換わります。1分たってもリモートライン 上にノットレディ状態が残っている場合は、このエラーメッセージが生成 されます。

考えられる原因

- リモートラインに接続されたいず ノットレディ状態になっている機器 れかの機器がノットレディ状態に が正しく設置され、分析に合わせて なっている。
 リモートケーブルの故障。
 リモートケーブルの故障。
 リモートケーブルを交換します。
- 3 ノットレディ状態になっている機
 その機器が故障していないか確認し
 器の部品の故障。
 ます(機器の付属書類を参照してく
 ださい)。



Shutdown

Error ID: 0063

シャットダウン

外部機器がリモートライン上にシャットダウンシグナルを生成しました。 モジュールは、リモート入力コネクタ上でステータスシグナルを常にモニ タしています。リモートコネクタのピン 4 に LOW シグナル入力がある と、このエラーメッセージが生成されます。

考えられる原因

- 1 システムへの CAN 接続により、別 外部機器内のリークを処理してから、 のモジュール内でリークが検出さ モジュールを再起動します。 れた。
- 2 システムへのリモート接続によ 外部機器内のリークを処理してから、 り、外部機器内でリークが検出さ モジュールを再起動します。 れた。
- 3 システムへのリモート接続によ 外部機器がシャットダウン状態に り、外部機器でシャットダウンが なっていないか確認します。 発生した。
Lost CAN Partner

Error ID: 0071

CAN 通信消失

分析中に、システム内の1台以上のモジュールの間で内部同期または通信 に失敗しました。

システムプロセッサは、システムコンフィグレーションを常にモニタリン グしています。1 台以上のモジュールとシステムの接続が認識されなくな ると、このエラーメッセージが生成されます。

考えられる原因

対策

1 CAN ケーブルの断線。

- すべての CAN ケーブルが正しく接続されていることを確認します。
- すべての CAN ケーブルが正しく設置されていることを確認します。
- 2 CAN ケーブルの不具合。 CAN ケーブルを交換します。
- 他のモジュールのメインボードのシステムをオフにします。システム 故障。
 を再起動して、システムが認識しないモジュールを確認します。



Timeout

Error ID: 0062

タイムアウト

タイムアウト値を超えました。

考えられる原因

- 1 分析が正常終了した後、要求通り ジュールをオフにしました。
- 2 シーケンスまたはマルチ注入測定 間、ノットレディ状態が続いた。

対策

ログブックを確認して、ノットレ にタイムアウト機能によってモ ディ状態が発生していないか、その 原因は何かを調べます。必要に応じ て、分析を再開してください。

ログブックを確認して、ノットレ 中に、タイムアウト値より長い時 ディ状態が発生していないか、その 原因は何かを調べます。必要に応じ て、分析を再開してください。

エラー情報 10 ポンプのエラー メッセージ

ポンプのエラー メッセージ

これらのエラーはポンプに固有のものです。

Encoder Missing

Error ID: 2046, 2050, 2510

エンコーダが見つからない

ポンプのポンプモータの光学エンコーダが見つからないか、故障していま す。

プロセッサは、ポンプエンコーダコネクタがあるかを 2 s ごとに確認しま す。プロセッサでコネクタが検出されないと、このエラーメッセージが生 成されます。

考えられる原因

- 1 ポンプエンコーダコネクタの不 Agilent Technologies に連絡してく ださい。 良、または接続されていない。
- 2 ポンプドライブアセンブリの不具 Agilent Technologies に連絡してく ださい。 合。

10 エラー情報 ポンプのエラー メッセージ

Index Adjustment

Error ID: 2204, 2214

インデックス調整

モジュールのエンコーダインデックス位置がズレています。

初期化中に、第1ピストンは機械的停止位置まで移動します。機械的停止 位置に到達後、エンコーダインデックスが到達するまで、ピストンは逆方 向に動きます。インデックス位置に到達する時間が長すぎると、このエ ラーメッセージが生成されます。

考えられる原因

対策

 ドライブの動きが不規則か、固着 している。
 ポンプヘッドを取り外し、シール、 ピストン、内部部品に摩耗、汚染、 損傷の兆候がないか検査します。必要に応じて、部品を交換します。
 ポンプドライブアセンブリの不具 合。
 Agilent Technologies に連絡してく ださい。

エラー情報 10 ポンプのエラー メッセージ

Index Limit

Error ID: 2203, 2213

インデックスリミット

ピストンがエンコーダインデックス位置に到達する所要時間が短すぎまし た (ポンプ)。

初期化中に、第1 ピストンは機械的停止位置まで移動します。機械的停止 位置に到達後、エンコーダインデックスが到達するまで、ピストンは逆方 向に動きます。インデックス位置に到達するのが早過ぎると、このエラー メッセージが生成されます。

考えられる原因

- 1 ドライブの動きが不規則か、固着 ポンプヘッドを取り外し、シール、 ピストン、内部部品に摩耗、汚染、 している。 損傷の兆候がないか検査します。必 要に応じて、部品を交換します。
- 2 ポンプドライブアセンブリの不具 Agilent Technologies に連絡してく ださい。 合。



Index Missing

Error ID: 2205, 2215, 2505

インデックスが見つからない

初期化中に、モジュールのエンコーダインデック位置が見つかりませんで した。

初期化中に、第1ピストンは機械的停止位置まで移動します。機械的停止 位置に到達後、エンコーダインデックスが到達するまで、ピストンは逆方 向に動きます。定義された時間内にインデックス位置が認識されないと、 このエラーメッセージが生成されます。

考えられる原因

- エンコーダケーブルの断線または Agilent Technologies に連絡してく 不具合。
- ポンプドライブアセンブリの不具 Agilent Technologies に連絡してく 合。

Initialization Failed

Error ID: 2207, 2217

初期化失敗

モジュールが、最大タイムウィンドウ内での初期化完了に失敗しまた。

ポンプ初期化サイクル全体に対して、最大時間が割り当てられます。初期 化完了前にこの時間を過ぎると、このエラーメッセージが生成されます。

ださい。

考えられる原因対策1 ポンプドライブアセンブリの不具
合。Agilent Technologies に連絡してく
ださい。2 メインボードの故障。Agilent Technologies に連絡してく

Missing Pressure Reading

Error ID: 2054

圧力測定値が読み取れない

ポンプの AD コンバータにより読み取られる圧力測定値が見つかりません。 AD コンバータは、ダンパーからの圧力シグナルを 1 ms ごとに読み取りま す。測定値が 10 s 以上不明の場合、エラーメッセージが生成されます。

考えられる原因

1 ダンパーの断線。 Agilent Technologies に連絡してください。

- **2** ダンパーの不具合。 Agilent Technologies に連絡してください。
- **3** メインボードの故障。 Agilent Technologies に連絡してください。

10 エラー情報 ポンプのエラー メッセージ

Motor-Drive Power

Error ID: 2041, 2042

モータドライブ電力

ポンプモータの消費電力が上限値を超過しました。

通常、流路の詰まりはダンパーの圧力センサーで検出され、圧力上限値を 超えるとポンプを電源を切ることになります。ダンパーの前で詰まりが発 生した、すなわち、圧力センサが圧力の上昇を検出できない場合は、モ ジュールは送液を続行します。圧力が上昇するにつれて、ポンプドライブ の消費電力はますます増えます。電流が上限値に達すると、モジュールの 電源が切れ、このエラーメッセージが生成されます。

考えられる原因

対策

ダンパーの前の流路が詰まってい ます。	ポンプヘッドとダンパーインレット 間のキャピラリとフリットが詰まっ ていないか確認します。
アウトレットバルブが詰まってい る。	アウトレットバルブを交換します。
ポンプドライブアセンブリの摩擦 が大きい (部分的な機械的障害物)。	ポンプヘッドアセンブリを取り外し ます。ポンプヘッドアセンブリまた はポンプドライブアセンブリの機械 的動作が妨害されていないことを確 認します。
ポンプドライブアセンブリの不具 合。	Agilent Technologies に連絡してく ださい。
	ダンパーの前の流路が詰まってい ます。 アウトレットバルブが詰まってい る。 ポンプドライブアセンブリの摩擦 が大きい (部分的な機械的障害物)。 ポンプドライブアセンブリの不具 合。

5 メインボードの故障。 Agilent Technologies に連絡してく ださい。

エラー情報 10 ポンプのエラー メッセージ

Pressure Above Upper Limit

Error ID: 2014, 2500

圧力が上限値を超過

システム圧力が圧力上限値を超過しました。

考えられる原因

対策

1	圧力上限値設定が低すぎます。	圧力上限が分析に適した値に設定さ れているか確認します。
2	流路に詰まりがある(<i>ダンパー</i> の 後)。	流路の詰まりを確認します。特に詰 まりやすい部分は次の部品です。イ ンラインフィルタフリット、ニード ル(オートサンプラ)、シートキャ ピラリ(オートサンプラ)、サンプ ルループ(オートサンプラ)、カラ ムフリット、細いキャピラリの使用

3 ダンパーの不具合。 Agilent Technologies に連絡してく ださい。

(内径 50 µm など)。

4 メインボードの故障。Agilent Technologies に連絡してく ださい。



Pressure Below Lower Limit

Error ID: 2015, 2501

圧力が下限値より低い

システム圧力が圧力下限値を下回りました。

考えられる原因

対策

1 圧力下限値設定が高すぎます。 圧力下限値が

2 リークがあります。

- 圧力下限値が分析に適した値に設定 されていることを確認します。
- ポンプヘッド、キャピラリ、 フィッティングにリークの兆候が ないかを検査します。
- モジュールをパージします。圧力 テストを実行して、シールまたは 他のモジュール部品に不具合がな いか確認します。
- **3** メインボードの故障。 Agilent Technologies に連絡してく ださい。

エラー情報 10 ポンプのエラー メッセージ

Pressure Signal Missing

Error ID: 2016

圧カシグナルが読み取れません

ダンパーの圧力シグナルが読み取れません。

ダンパーの圧力シグナルは、指定電圧範囲内にある必要があります。圧力 シグナルが失われた場合、プロセッサは、ダンパーコネクタを通じて、約 -120 mV の電圧を検出します。

考	えられる原因	対策		
1	ダンパーの断線。	Agilent Technol	logies に連絡してください	•
2	ダンパーの不具合。	Agilent Technol	logies に連絡してください	۰ _°

Pump Configuration

Error ID: 2060

ポンプ コンフィグレーション

電源を入れた時に、グラジエントポンプが新しいポンプ コンフィグレー ションを認識しました。

グラジエントポンプは、工場でコンフィグレーションを割り当てられます。 グラジエントバルブ切断してグラジェントポンプを再起動すると、エラー メッセージを発します。しかし、ポンプはこのコンフィグレーションでア イソクラティックポンプとして機能します。電源を入れるたびにエラー メッセージが再表示されます。

考えられる原因

対策

 グラジエントバルブが接続されて がない。
 グラジエントバルブを再び接続しま す。 **10 エラー情報** ポンプのエラー メッセージ

Pump Head Missing

Error ID: 2202, 2212

ポンプヘッドが見つからない

ポンプのポンプヘッドの末端位置が見つかりませんでした。

ポンプを再起動すると、調整ドライブが機械的末端位置まで前進します。 通常は 20 s 以内に末端位置に到達します。これは、モータ電流の上昇で 示されます。20 s 以内に末端位置が見つからない場合、このエラーメッ セージが生成されます。

考えられる原因

対策

- ポンプヘッドが正しく取り付けら れていない(ネジが固定されてい ないか、ポンプヘッドがしっかり 固定されていない)。
 ポンプヘッドを正しく取り付けます。
 ポンプヘッドを正しく取り付けます。
- 2 ピストンの破損。

ピストンを交換します。

Restart Without Cover

Error ID: 2502

カバーなし再スタート

上部カバーと発泡材が開いた状態でモジュールが再起動されました。

上部発泡材が定位置にくると、メインボード上のセンサーによって検出さ れます。発泡材を取り外した状態でモジュールを再起動すると、モジュー ルは 30 秒以内に電源が切れ、このエラーメッセージが生成されます。

考えられる原因

対策

1 上部カバーおよび発泡材を取り外した Agilent のサービス担当者に連 状態でモジュールを起動しました。 絡してください。

Servo Restart Failed

Error ID: 2201, 2211

サーボ再起動の失敗

モジュールのポンプモータが、再起動時の正しい位置まで動くことができませんでした。

モジュールの電源を入れると、最初の段階で可変磁気抵抗モータの C 相が オンになります。このとき、ロータは C 位置の 1 つに移動する必要があ ります。サーボが整流子とのフェーズシーケンス処理をコントロールでき るようにするためには、ロータが C 位置にある必要があります。ロータが 動くことができないか、C 位置に到達しない場合は、このエラーメッセー ジが生成されます。

考えられる原因

対策

1	ケーブルの断線または不具合。	Agilent Technologies に連絡してく ださい。
2	モジュールの機械的動作の妨害。	ポンプヘッドアセンブリを取り外し ます。ポンプヘッドアセンブリまた はポンプドライブアセンブリの機械 的動作が妨害されていないことを確 認します。
3	ポンプドライブアセンブリの不具 合。	Agilent Technologies に連絡してく ださい。

4 メインボードの故障。Agilent Technologies に連絡してく ださい。



Stroke Length

Error ID: 2206, 2216

ストローク長

ピストンの下限位置と機械的上端位置の距離がリミット値から外れていま す (ポンプ)。

初期化中に、モジュールはドライブ電流をモニタリングします。ピストン が予想よりも早く上の機械的末端位置に達した場合、モジュールはピスト ンを末端位置の先まで押し上げようとするため、モータ電流が上がります。 この電流が上昇すると、このエラーメッセージが生成されます。

考えられる原因

対策

 ポンプドライブアセンブリの不具 Agilent Technologies に連絡してく 合。
 ださい。

エラー情報 10 ポンプのエラー メッセージ

Temperature Limit Exceeded

Error ID: 2517

温度上限を超過

モータドライブ回路の1つの温度が高過ぎます。

プロセッサは、メインボード上のドライブ回路の温度を常にモニタリング しています。長時間にわたって消費電流が大きすぎると、ドライブ回路の 温度は上昇します。この温度が上限値を超えると、このエラーメッセージ が生成されます。

考えられる原因

1	ポンプドライブアセンブリの摩擦 が大きい (部分的な機械的障害物)。	ポンプヘッドアセンブリを取り外し ます。ポンプヘッドアセンブリまた はポンプドライブアセンブリの機械 的動作が妨害されていないことを確 認します。
2	ダンパーの前の流路が部分的に詰 まっています。	アウトレットバルブが詰まっていな いことを確認します。
3	ポンプドライブアセンブリの不具 合。	Agilent Technologies に連絡してく ださい。
4	メインボードの故障。	Agilent Technologies に連絡してく ださい。



Temperature Out of Range

Error ID: 2517

範囲外の温度

モータドライブ回路の温度センサ測定値が範囲外です。

ハイブリッドセンサから ADC に供給される値は $0.5 V \sim 4.3 V$ の範囲内 でなければなりません。値がこの範囲を外れると、このエラーメッセージ が生成されます。

考えられる原因

対策

1 メインボードの故障。 Agilent Technologies に連絡してください。

エラー情報 10

ポンプのエラー メッセージ

Valve Failed (MCGV)

Error ID: 2040

バルブ故障 (MCGV)

- バルブ 0 故障:バルブ A
- バルブ1 故障:バルブ B
- バルブ 2 故障:バルブ C
- バルブ 3 故障:バルブ D

複数のチャンネルグラジエントバルブの内の 1 つのバルブが、正しく切り 換えられなかつた。

プロセッサが、各切り換えサイクル前後のバルブ電圧をモニタリングして います。電圧が予想リミット値を外れると、このエラーメッセージが生成 されます。

考えられる原因

- グラジエントバルブが接続されて グラジエントバルブが正しく接続さいない。
 れているか確認します。
 (機器内部の)接続ケーブルが接 Agilent Technologies に連絡してく
- 2 (機器Piplo) 接続の ジルが後 Agrient Technologies に座格 じてく 続されていません。 ださい。
- **3** (機器内の)接続ケーブルの不具 Agilent Technologies に連絡してく 合。 ださい。
- 4 グラジエントバルブの不具合。 グラジエントバルブを交換します。

10 エラー情報 ポンプのエラー メッセージ

MCGV Fuse

Error ID: 2043

MCGV ヒューズ

バルブヒューズ 0: チャンネル A と B

バルブヒューズ 1: チャンネル C と D

クォータナリポンプのグラジエントバルブの消費電力が大きすぎたため、 その電子ヒューズが切れました。

考えられる原因

- グラジエントバルブの不具合。 エラーメッセージが再度表示された 場合、グラジエントバルブを交換し ます。

 接続ケーブル(正面パネルからメ インボードまで)の不具合。
 Agilent のサービス担当者に連絡し てください。
- **3** メインボードの故障。 Agilent Technologies に連絡してく ださい。

エラー情報 10

ポンプのエラー メッセージ

Wait Timeout

Error ID: 2053

タイムアウト待ち

診断モードまたは特殊なアプリケーションで特定のテストを実行する場合、 ポンプは、ピストンが特定の位置に到達するまで待機するか、特定の圧力 または流量に到達するまで待機する必要があります。それぞれの動作また は状態はタイムアウト期間内に完了する必要があります。そうでなかった 場合には、エラーメッセージが発せられます。

完了しないと、このエラーメッセージが生成されます。

圧力が到達しなかった。

ポンプチャンネル A が送液段階に到達しなかった。

ポンプチャンネル B が送液段階に到達しなかった。

ポンプチャンネル A が吸引段階に到達しなかった。

ポンプチャンネル B が吸引段階に到達しなかった。

指定時間内に溶媒容量が送液されなかった。

考えられる原因

- 1 テストの開始後、流量が変化し 使用中の特殊なアプリケーションに 適した動作条件か確認します。 た。
- 2 ポンプドライブアセンブリの不具 Agilent のサービス担当者に連絡し てください。 合。



Solvent Zero Counter

Error ID: 2055, 2524

溶媒ゼロカウンタ

ポンプファームウェアバージョン A. 02. 32 以降により、データシステムで 溶媒ボトル補充の設定が可能になります。機能が設定されている場合、そ れに応じてボトル中の容量レベルが指定値より下だと、エラーメッセージ が表示されます。

考えられる原因

- ボトルの容量が指定容量を下回り ボトルを再補充して、溶媒カウンタ ました。
 をリセットします。
- 設定が正しくありません。
 リミット値が正しく設定されている か確認してください。

オート サンプラのエラー メッセージ

これらのエラーはオート サンプラに固有のものです。

Arm Movement Failed

Error ID: 4002

アーム移動失敗

トランスポートアセンブリが、いずれかの軸の動作を完了できませんでした。

プロセッサでは、個々の軸方向での動作が正常に完了するまでのタイム ウィンドウを定義しています。トランスポートアセンブリの動作とポジ ションは、ステッピングモーター上のエンコーダによってモニタリングさ れます。プロセッサがこのエンコーダから正確なポジション情報をタイム ウィンドウ内で受信しないと、このエラーメッセージが生成されます。

それぞれの軸については、『109ページ図 37』の図を参照してください。

- ・ [アーム移動失敗 0]: X 軸
 - [アーム移動失敗 1]: Z 軸
 - [アーム移動失敗 2]:シータ(グリッパの回転)
 - [アーム移動失敗 3]: グリッパ (グリッパフィンガーの開閉)

考えられる原因	対策
1 機械的に妨害を受けている。	トランスポートアセンブリの動作が 阻害されていないか、確認します。
 トランスポートアセンブリの摩擦 が大きい。 	Agilent のサービス担当者に連絡し てください。
3 モータアセンブリの故障。	Agilent のサービス担当者に連絡し てください。
4 サンプル トランスポート アセン ブリのフレックスボードの故障。	Agilent のサービス担当者に連絡し てください。
5 メインボードの故障。	Agilent のサービス担当者に連絡し てください。

Initialization Failed

Error ID: 4020

初期化失敗

オートサンプラは初期化を正常に完了できませんでした。

オートサンプラの初期化作業は、ニードルアームとトランスポートアセン ブリを、定義済みのシーケンスで、それらのホーム位置に移動します。初 期化中、プロセッサは、ポジションセンサとモーターエンコーダの動作が 正しいかモニタリングします。動作が正常に終了しない、または検出され ないと、このエラーメッセージが生成されます。

考えられる原因

1	機械的に妨害を受けている。	トランスポートアセンブリの動作が 阻害されていないか、確認します。
2	サンプリングユニットのフレック スボードの故障。	Agilent のサービス担当者に連絡し てください。
3	トランスポートアセンブリのフ レックスボードの故障。	Agilent のサービス担当者に連絡し てください。
4	サンプリングユニットのモータの 故障。	Agilent のサービス担当者に連絡し てください。
5	メインボードの故障。	Agilent のサービス担当者に連絡し てください。

Initialization with Vial

Error ID: 4028

バイアル初期化

オートサンプラが、グリッパ内にまだバイアルがある間に初期化しようとしました。

初期化中、オートサンプラは、モーターエンコーダのモニタ中にグリッパ フィンガーを開閉することによって、グリッパを正常に操作できるかどう かをチェックします。初期化が開始されたとき、グリッパ内にまだバイア ルがある場合には、グリッパフィンガーは閉じず、エラーメッセージが生 成されます。

考えられる原因

対策

 グリッパ内にまだバイアルがあ る。
 ユーザーインターフェースの リリー
 スバイアル機能を使用して、バイア ルを取り除きます。オートサンプラ

ルを取り除きます。オートサンプラを再初期化します。

Invalid Vial Position

Error ID: 4042

無効なバイアルポジション

メソッドまたはシーケンス内で設定されているバイアルポジションが存在 しません。

どのサンプルトレイが取り付けられているかは、トランスポートアセンブ リのフレックスボード上にある反射センサーにより自動的にチェックされ ます。バイアルのポジションが現在のサンプルトレイのコンフィグレー ション内に存在しないと、このエラーメッセージが生成されます。

考えられる原因

対策

- 間違ったトレイがインストールさ 正しいトレイを取り付けるか、メ れている。
 ソッドまたはシーケンスを変更しま す。
- 2 メソッドまたはシーケンス内で設 取り付けれらたトレイがメソッド設定されているバイアルポジション 定と一致しているかを確認します。 が間違っている。
- 3 トレイの認識の誤り(サンプルト サンプルトレイの背面にあるコーレイが汚れている、またはトランディング表面が汚れていないかを確スポートアセンブリのフレックス認します。問題が解決しない場合は、ドランスポートアセンブリを交換し

てください。

Metering Home Failed

Error ID: 4054, 4704

計量ピストンのホームポジションへの移動失敗

計量ピストンが、ホームポジションに戻りませんでした。

計量ピストンのホームポジションは、サンプリングユニットのフレックス ボード上にあるホームポジションセンサーでモニタリングされます。プラ ンジャがホームポジションに戻らなかったり、センサーがプランジャのポ ジションを認識できないと、このエラーメッセージが生成されます。

考	えられる原因	対策
1	センサーが汚れているか、故障し ている。	Agilent のサービス担当者に連絡し てください。
2	ピストンの破損。	メタリングピストンとシールを交換 します。
3	計量ドライブのモータの故障。	Agilent のサービス担当者に連絡し てください。
4	メインボードの故障。	Agilent のサービス担当者に連絡し てください。

Missing Vial

Error ID: 4019, 4034, 4541, 4706

バイアルなし

メソッドまたはシーケンス内で設定されているポジションにバイアルが見 つかりませんでした。

グリッパアームがサンプルトレイからバイアルを取り上げるとき、プロ セッサは、グリッパモーターのエンコーダをモニタします。バイアルが存 在する場合には、グリッパフィンガーを閉じる動作は、バイアルによって 制限されます。一方、バイアルが存在しない場合には、グリッパフィン ガーの閉じる動作が大きくなります。これはプロセッサによって検出され (エンコーダの位置)、エラーメッセージが生成されます。

考えられる原因

- 1 メソッドまたはシーケンス内で設 サンプルバイアルを正しいポジショ ンに置くか、あるいはメソッドまた 定されているポジションにバイア はシーケンスを変更します。 ルがない。
- 2 グリッパのポジション調節不良。 グリッパのポジションを調節します。 (アライメント不良)
- 3 グリッパアセンブリの不良 (グ グリッパアセンブリを交換します。 リッパの指またはベルトの不良)。
- **4** トランスポートアセンブリのフ Agilent のサービス担当者に連絡し てください。 レックスボードの故障。

Missing Wash Vial

Error ID: 4035, 4542, 4707

洗浄用バイアルが見つからない

メソッドでプログラムされた洗浄用バイアルが見つかりません。

グリッパアームがサンプルトレイからバイアルを取り上げるとき、プロ セッサは、グリッパモーターのエンコーダをモニタします。バイアルが存 在する場合には、グリッパフィンガーを閉じる動作は、バイアルによって 制限されます。一方、バイアルが存在しない場合には、グリッパフィン ガーの閉じる動作が大きくなります。これはプロセッサによって検出され (エンコーダの位置)、エラーメッセージが生成されます。

考えられる原因

対策

 メソッドで定義された位置に洗浄 洗浄用バイアルを正しいポジション バイアルがない。
 に置くか、メソッドを変更します。

Motor Temperature

Error ID: 4027, 4040, 4261, 4451

モーター温度

トランスポートアセンブリのモーターに過度の電流が流れたため、その モーターが高温になっています。プロセッサは、そのモーターが破損しな いように、そのモーターの電源を切りました。

個々のモーターについては 『109 ページ 図 37』の図を参照してください。

- [モーター温度 0]: X 軸モーター
 - 「**モーター**温度 11: Z 軸モーター
 - [モーター温度 2]:シータ軸(グリッパの回転)モーター
 - [モーター温度 3]: グリッパモーター (グリッパフィンガー用 モーター)

プロセッサは、各モーターに流れる電流とモーターに電流が流れた時間を モニタリングします。一連のモーターに流れる電流は、各モーターの負荷 (摩擦、部品の大きさなど)によって決まります。電流が高すぎたり、電 流がモーターに流れる時間が長すぎた場合に、このエラーメッセージが生 成されます。

考えられる原因

対策

1	機械的に妨害を受けている。	トランスポートアセンブリの動作が阻 害されていないか、確認します。
2	トランスポートアセンブリの摩 擦が大きい。	Agilent のサービス担当者に連絡して ください。
3	モータのベルトの張りが強すぎ る。	電源スイッチでオートサンプラをオフ にします。最低 10 min 待機してから、 スイッチをもう一度オンにします。
4	モータの故障。	Agilent のサービス担当者に連絡して ください。
Б	トランスポートアセンブルのフ	Agilant のサービス担当者に演練して

5 トランスポートアセンブリのフ Agilent のサービス担当者に連絡して レックスボードの故障。 ください。

Needle Down Failed

Error ID: 4018

ニードルダウンの失敗

ニードルアームが、ニードルシートに降りることができませんでした。

ニードルアームの下方位置は、サンプリングユニットのフレックスボード 上の位置センサーによって、モニタされています。センサーは、ニードル シート位置への移動が完了すると、それを検出します。ニードルが終了位 置に達しなかった、またはセンサーがニードルアームの移動を検出できな かった場合には、エラーメッセージが生成されます。

考えられる原因

対策

- ニードルが正しくインストールさ 正しいニードルタイプを使用していれていない。またはニードルタイ て、正しくインストールされているプが間違っている(長すぎる)。ことを確かめます。
- 位置センサーの不良または汚れ。 Agilent のサービス担当者に連絡してください。

てください。

Agilent のサービス担当者に連絡し

3 モータの故障。

- 4 スピンドルアセンブリが引っか Agilent のサービス担当者に連絡しかっている。
- 5 メインボードの故障。 Agilent のサービス担当者に連絡し てください。

Needle Up Failed

Error ID: 4017

ニードルアップの失敗

ニードルアームが、シートから、またはバイアルから上部位置に正常に移動しませんでした。

ニードルアームの上部位置は、サンプリングユニットのフレックスボード 上の位置センサーによって、モニタされています。センサーは、ニードル の上部位置への移動が完了すると、それを検出します。ニードルが終了位 置に達しなかった、またはセンサーがニードルアームの移動を検出できな かった場合には、エラーメッセージが生成されます。

考えられる原因

1	位置センサーの不良または汚れ。	Agilent のサービス担当者に連絡し てください。
2	モータの故障。	Agilent のサービス担当者に連絡し てください。
3	スピンドルアセンブリが引っか かっている。	Agilent のサービス担当者に連絡し てください。
4	メインボードの故障。	Agilent のサービス担当者に連絡し てください。

Safety Flap Missing

Error ID: 4032

安全フラップ不良

安全フラップが検出されませんでした。

サンプルを注入するためにニードルがニードルシートに降りる前に、安全 フラップは定位置にロックされます。次に、グリッパは安全フラップを ニードルから除去する動作で安全フラップをチェックします。グリッパが 安全フラップの位置を越えて移動することができた (安全フラップが位置 にない)場合には、エラーメッセージが生成されます。

考えられる原因

対策

安全フラップがない、または壊れ Agilent のサービス担当者に連絡している。
 てください。

Valve to Bypass Failed

Error ID: 4014, 4701

バイパスへのバルブ切り替え失敗

インジェクションバルブをバイパスポジションに切り替えることができま せんでした。

インジェクションバルブの切り替えは、バルブアセンブリ上にある 2 つの マイクロスイッチによってモニタリングされます。これらのスイッチで、 インジェクションバルブの動作が正常に完了したかを検出します。イン ジェクションバルブがバイパスポジションに到達できないか、あるいはマ イクロスイッチが閉じないと、このエラーメッセージが生成されます。

考えられる原因

1	インジェクションバルブの故障。	Agilent のサービス担当者に連絡し てください。
2	メインボードの故障。	Agilent のサービス担当者に連絡し てください。

Valve to Mainpass Failed

Error ID: 4015

メインパスへのバルブ切り替え失敗

インジェクションバルブをメインパスポジションに切り替えることができ ませんでした。

インジェクションバルブの切り替えは、バルブアセンブリ上にある 2 つの マイクロスイッチによってモニタリングされます。これらのスイッチで、 インジェクションバルブの動作が正常に完了したかを検出します。イン ジェクションバルブがメインパスポジションに到達できないか、あるいは マイクロスイッチが閉じないと、このエラーメッセージが生成されます。

考えられる原因

1	インジェクションバルブの故障。	Agilent のサービス担当者に連絡し てください。
2	メインボードの故障。	Agilent のサービス担当者に連絡し てください。

Vial in Gripper

Error ID: 4033

グリッパ内にバイアルあり

グリッパアームが、まだグリッパ内にバイアルがある間に移動しようとし ました。

サンプリング動作の特定のステージでは、グリッパがバイアルを保持する ことはできません。オートサンプラは、モーターエンコーダのモニタ中に、 グリッパフィンガーを開閉することによって、サンプルのバイアルがグ リッパ内にあるかどうかをチェックします。グリッパフィンガーが閉じる ことができなかった場合には、エラーメッセージが生成されます。

考えられる原因

対策

1 グリッパ内にまだバイアルがあ ユーザーインターフェースの リリー **スバイアル**機能を使用して、バイア る。 ルを取り除きます。オートサンプラ

を再初期化します。

10 エラー情報

一般検出器エラー メッセージ

一般検出器エラー メッセージ

これらのエラーは、VWD および DAD の両方の検出器タイプに固有のものです。

Heater at fan assembly failed

Error ID: 1073

ファンアセンブリのヒーター故障

重水素ランプまたはタングステンランプ(DAD のみ)の電源をオン / オフ を行うごとに、ヒーターの自己診断が行われます。テストが失敗すると、 エラーイベントが作成されます。結果として、温度コントロールはオフに なります。

考えられる原因

対策

コネクタまたはケーブルの不良。 Agilent のサービス担当者に連絡してください。
 ヒータが故障しています。 Agilent のサービス担当者に連絡してください。
エラー情報 10 た山岡エニー メッセージ

一般検出器エラー メッセージ

Heater Power At Limit

Error ID: 1074

ヒーター出力限界

ヒーターの出力が上限値または下限値に到達しました。このイベントは、 運転あたり 1 回のみ送信されます。パラメータにより、どのリミット値を ヒットしたかが判ります。

0 は出力上限値ヒットを意味します(過剰な周囲温度降下)。

1 は出力下限値ヒットを意味します(過剰な周囲温度上昇)。

考えられる原因

対策

1 周囲温度の過度な変化 温度コントロールが平衡化するまで 待機します。

Illegal temperature value from sensor at fan assembly

Error ID: 1071

ファンアセンブリのセンサからの不正な温度値

温度センサが許容範囲外の値を送信しました。このイベントのパラメータ は、測定温度と 1/100 ℃で一致します。結果として、温度コントロールは オフになります。

考えられる原因

- センサーが汚れているか、故障し Agilent のサービス担当者に連絡している。
 てください。
- 検出器が不正な周囲条件に曝され 周囲条件が許容範囲内であることを ています。
 確認します。

Illegal Temperature Value from Sensor at Air Inlet

Error ID: 1072

空気吸入口のセンサからの不正な温度値

温度センサ(検出器メインボードに設置)が、許容範囲外の値を送信しま した。このイベントのパラメータは、測定温度と 1/100 ℃で一致します。 結果として、温度コントロールはオフになります。

考えられる原因

ています。

1	温度センサが故障しています。	Agilent のサービス担当者に連絡し てください。
2	検出器が不正な周囲条件に曝され ています。	周囲条件が許容範囲内であることを 確認します。

エラー情報 10

一般検出器エラー メッセージ

UV Lamp Current

Error ID: 7450

UV ランプ電流

UV ランプ電流がありません。

運転中、プロセッサはランプに取り込まれるアノード電流を継続的にモニ タリングします。アノード電流が下限値を下回ると、このエラーメッセー ジを発します。

考えられる原因

- ランプが外れている。
 UW ランプコネクタがしっかりと固定 されているか確認します。
- **2** UV ランプの故障またはアジレン UV ランプを交換します。 ト以外のランプ。
- 6 検出器メインボードの故障。
 7 ジレントのサービス担当者にご連絡ください。
- 4 電源の故障。 アジレントのサービス担当者にご連絡ください。

10 エラー情報 一般検出器エラー メッセージ

UV Lamp Voltage

Error ID: 7451

UV ランプ電圧

UV ランプアノード電圧がありません。

運転中、プロセッサはランプのアノード電圧を継続的にモニターします。 アノード電圧が下限値を下回ると、このエラーメッセージを発します。

考えられる原因

- UV ランプの故障またはアジレン UV ランプを交換します。
 ト以外のランプ。
- 2 検出器メインボードの故障。 おください。

 3 電源の故障。
 アジレントのサービス担当者にご連
 アジレントのサービス担当者にご連
 - アジレントのサービス担当者にご連 絡ください。

エラー情報 10 た出業エラーメッセージ

VWD 検出器エラーメッセージ

VWD 検出器エラーメッセージ

これらのエラーは可変波長検出器に固有のものです。

ADC Hardware Error

Error ID: 7830, 7831

ADC ハードウェアエラー

A/D コンバータハードウェアが故障しています。

考えられる原因

対策

1 A/D コンバータハードウェアが故 Agilent のサービス担当者に連絡し 障しています。 てください。

10 エラー情報 WWD 検出器エラーメッセージ

Wavelength calibration setting failed

Error ID: 7310

波長キャリブレーション設定の失敗

波長キャリブレーション中に最大強度が検出されませんでした。

キャリブレーション 0 失敗:	ゼロオーダキャリブレーションに失 敗しました。
キャリブレーション 1 失敗:	656 nm キャリブレーションに失敗し ました。

考えられる原因

ズレ / 故障。

- 1 ランプが消灯している。 ランプの電源を入れます。
- 2 フローセルの設置が間違ってい フローセルが正しく設置されている ことを確認します。 る。
- 3 フローセルの汚れや気泡がある。 フローセルウィンドウのクリーニン グ / 交換または気泡の除去を行いま す。
- 4 強度が低すぎる。 ランプを交換します。
- 5 現在のステップ値が最大値から大 ・ キャリブレーションを繰り返しま きく離れている。 す。
 - Agilent のサービス担当者に連絡 してください。
- 6 グレーティングアセンブリの位置 Agilent のサービス担当者に連絡し てください。
- 7 メインボードの故障。 Agilent のサービス担当者に連絡し てください。

Cutoff filter doesn't decrease the light intensity at 250 nm

Error ID: 7813

250 nm でカットオフフィルタによる光強度の減少なし

ランプ点火後の自動フィルタチェックに失敗しました。ランプ点灯後、検 出器はカットオフフィルタを光路の中に移動させます。フィルタが正しく 機能していると、ランプ強度の減少が見られます。予想される強度減少が 検出されないと、このエラーメッセージが生成されます。

考えられる原因		対策	
1	モーターが接続されていません。	Agilent のサービス担当者に連絡し てください。	
2	モーターの故障	Agilent のサービス担当者に連絡し てください。	
3	グレーティングまたはフィルタの 故障 / 消失	Agilent のサービス担当者に連絡し てください。	
4	ケーブル / コネクタの不具合	Agilent のサービス担当者に連絡し てください。	

10 エラー情報 VWD 検出器エラーメッセージ

Filter Missing

Error ID: 7816

フィルタがありません

フィルタ モータが検出されません。

考えられる原因

- フィルタモータが接続されていま Agilent のサービス担当者に連絡し せん。
- **2** ケーブル / コネクタの不具合。 Agilent のサービス担当者に連絡してください。

エラー情報 10 WWD 検出器エラーメッセージ

Grating or Filter Motor Errors

Error ID: Grating: 7800, 7801, 7802, 7803, 7804, 7805, 7806, 7808, 7809; Filter: 7810, 7811, 7812, 7813, 7814, 7815, 7816

グレーティング / フィルタモーターエラー

モーターテストに失敗しました。

- テスト 0 失敗: フィルタモーター
- グレーティングモーター テスト1 失敗:

モーターテスト中、検出器はエンドポジションセンサをモニタリングしな がら、モーターをエンドポジションに移動させます。エンドポジションが 検出されないと、このエラーメッセージが生成されます。

考えられる原因		対策	
1	モーターが接続されていません。	Agilent のサービス担当者に連絡し てください。	
2	モーターの故障	Agilent のサービス担当者に連絡し てください。	
3	グレーティングまたはフィルタの 故障 / 消失	Agilent のサービス担当者に連絡し てください。	
4	ケーブル / コネクタの不具合	Agilent のサービス担当者に連絡し てください。	

10 エラー情報 VWD 検出器エラーメッセージ

Grating Missing

Error ID: 7819

グレーティングがありません

グレーティングモータが検出されません。

考えられる原因

- グレーティングモータが接続され Agilent のサービス担当者に連絡していません。
- **2** ケーブル / コネクタの不具合。 Agilent のサービス担当者に連絡してください。

エラー情報 10 た山際エニーメッカージ

WWD 検出器エラーメッセージ

No heater current

Error ID: 7453

ヒーター電流なし

検出器のランプヒーター電流が流れていません。ランプ点火中、プロセッ サはヒーター電流をモニタリングします。電流が1 以内に下限値を超え ないと、このエラーメッセージが生成されます。

考えられる原因

1	ランプが外れている。	ランプが接続されているかを確認し ます。
2	上部発泡材が所定の位置にない状 態で、点火が開始された。	Agilent のサービス担当者に連絡し てください。
3	メインボードの故障。	Agilent のサービス担当者に連絡し てください。
4	故障か Agilent 製以外のランプ	ランプを交換します。
5	電源の故障。	Agilent のサービス担当者に連絡し てください。

Wavelength holmium check failed

Error ID: 7318

波長ホルミウムチェックの失敗

検出器のホルミウムオキサイドテストに失敗しました。ホルミウムオキサ イドテスト中、検出器はホルミウムフィルタを光路の中に移動させ、ホル ミウムオキサイドフィルタの測定吸光度の最大値と予想最大値を比較しま す。測定最大値がリミット値を外れると、このエラーメッセージが生成さ れます。

考えられる原因

1 グレーティングアセンブリの位置 ズレ / 故障。

- フローセルが正しく取り付けられていること、および汚染物質がないこと(セル窓、緩衝液など)を確認します。
- フィルタモーターテストを実行 し、フィルタモーターアセンブリ が故障していないかを確認しま す。故障している場合は、 Agilentのサービス担当者に連絡 してください。
- グレーティングモーターテストを 実行し、グレーティングアセンブ リが故障していないかを確認しま す。故障している場合は、 Agilentのサービス担当者に連絡 してください。

エラー情報 10 WWD 検出器エラーメッセージ

Ignition Failed

Error ID: 7452

点火不良

ランプが点火に失敗しました。点火サイクル中、プロセッサはランプ電流 をモニタリングします。ランプ電流が 2-5 s 以内に下限値を超えない と、このエラーメッセージが生成されます。

考えられる原因

1	ランプが外れている。	ランプが接続されているかを確認し ます。
2	故障か Agilent 製以外のランプ	ランプを交換します。
3	電源の故障。	Agilent のサービス担当者に連絡し てください。
4	メインボードの故障。	Agilent のサービス担当者に連絡し てください。



Wavelength test failed

Error ID: 7890

波長テスト失敗

ランプ点火後の自動波長チェックに失敗しました。ランプを点灯すると、 検出器はランプのウォームアップのために 1 min 待機します。その後、リ ファレンスダイオードによる重水素エミッションライン (656 nm) の チェックを行います。エミッションラインが 656 nm から 3 nm 以上離れ ていると、このエラーメッセージが生成されます。

考えられる原因

対策

正しくキャリブレーションされ 検出器のリキャリブレーションを行なった。
 います。

エラー情報 10
たりピュー イッセージ

DAD 検出器エラー メッセージ

DAD 検出器エラー メッセージ

これらのエラーはダイオードアレイ検出器に固有のものです。

可視ランプ電圧

可視ランプ電圧がありません。

運転中、プロセッサはランプ電圧を継続的にモニタリングします。電圧が 下限値を下回ると、このエラーメッセージが生成されます。

考えられる原因

対策

コネクタまたはケーブルの不良。 Agilent のサービス担当者に連絡してください。
 電源の故障。 Agilent のサービス担当者に連絡してください。



Diode Current Leakage

Error ID: 1041

ダイオード電流リーク

検出器の電源を入れると、プロセッサが各光学系ダイオードのリーク電流 を確認します。リーク電流が上限値を超えると、このエラーメッセージを 発します。

考えられる原因

1	PDA/ 光学ユニットの故障。	アジレントのサービス担当者にご連 絡ください。
2	コネクタまたはケーブルの不良。	アジレントのサービス担当者にご連 絡ください。

エラー情報 10 DAD 検出器エラー メッセージ

UV Ignition Failed

Error ID: 7452

UV 点灯失敗

UV ランプが点火に失敗しました。

点灯サイクル中、プロセッサが UV ランプ電流をモニターします。ランプ 電流が 2 ~ 5 秒以内に下限値を上回らないと、このエラーメッセージを 発します。

考えられる原因

対策

- 1 ランプが熱すぎます。高温ガス放 ランプを消灯し、冷却するまで 15 電ランプは、低温ランプほど簡単 分以上放置します。 に点火しないかもしれません。 2 ランプが外れている。
 - ランプが接続されているか確認しま す。
- **3** UV ランプの故障またはアジレン UV ランプを交換します。 ト以外のランプ。
- 4 検出器メインボードの故障。 アジレントのサービス担当者にご連 絡ください。 アジレントのサービス担当者にご連 5 電源の故障。

絡ください。

10 エラー情報

DAD 検出器エラー メッセージ

UV Heater Current

Error ID: 7453

UV ヒーター電流

UV ランプヒーター電流がありません。

UV ランプ点灯中、プロセッサはヒーター電流をモニターします。電流が 1 秒以内に下限値を上回らないと、このエラーメッセージを発します。

考えられる原因

1	ランプが外れている。	UV ランプが接続されているか確認し ます。
2	上部発泡材が所定の位置にない状 態で、点火が開始された。	アジレントのサービス担当者にご連 絡ください。
3	UV ランプの故障またはアジレン ト以外のランプ。	UV ランプを交換します。
4	検出器メインボードの故障。	アジレントのサービス担当者にご連 絡ください。
5	電源の故障。	アジレントのサービス担当者にご連 絡ください。

エラー情報 10 DAD 検出器エラー メッセージ

Calibration Values Invalid

Error ID: 1036

キャリブレーション値無効

分光計 ROM から読み取られたキャリブレーション値は無効です。

リキャリブレーション後、キャリブレーション値は ROM に保存されます。 プロセッサにより、キャリブレーションデータが有効かを定期的にチェッ クします。データが無効か、分光計 ROM から読み取ることができない場 合、エラーメッセージを発します。

考えられる原因 対策 1 コネクタまたはケーブルの不良。 アジレントのサービス担当者にご連絡ください。 2 PDA/光学ユニットの故障。 アジレントのサービス担当者にご連絡ください。



Holmium Oxide Test Failed

ホルミウムオキサイドテスト失敗

考えられる原因

1	ランプが消灯している。	ランプが点灯していることを確認し ます。
2	フローセルの故障または汚れ。	フローセルが正しく挿入されている こと、および汚染物質がないこを (セル窓、緩衝液など)を確認しま す。
3	フィルターアセンブリの故障。	Agilent のサービス担当者に連絡し てください。
4	アクロマートアセンブリの故障。	Agilent のサービス担当者に連絡し てください。
5	PDA/ 光学ユニットの故障。	Agilent のサービス担当者に連絡し てください。

エラー情報 10 DAD 検出器エラー メッセージ

Wavelength Recalibration Lost

Error ID: 1037

波長リキャリブレーション消失

検出器を正しく操作するために必要なキャリブレーション情報が失われました。

検出器のキャリブレーション中、キャリブレーション値は ROM に保存され ます。分光光度計 ROM にデータがないと、このエラーメッセージを発しま す。

考えられる原因		対策	
1	検出器が新品。	検出器のリキャリブレーションを行 います。	
2	検出器を修理した。	アジレントのサービス担当者にご連	

アジレントのサービス担当者にご連 絡ください。

10 エラー情報 DAD 検出器エラー メッセージ

DSP Not Running

DSP が実行されていません

光学ユニットとメインボードの間の通信に問題がある場合、このエラー メッセージが表示されます。

考えられる原因

- ランダムな通信エラー。
 電源スイッチで検出器の電源を切り、再び入れます。エラーが再発する場合、以下の操作を行います。
 Agilent のサービス担当者に連絡してください。
- 2 検出器メインボードの故障。 Agilent のサービス担当者に連絡してください。
 3 PDA/光学ユニットの故障。 Agilent のサービス担当者に連絡し
 - Agrient のリーヒス担当有に連絡し てください。

No Run Data Available In Device

デバイスに使用可能な分析データなし

非常にまれに、コンパクトフラッシュカードの容量が十分ではありません。 たとえば LAN 通信の遮断が長く、検出器が特別な設定を使用する場合(80 Hz のフルデータレートに加え、フルスペクトルとすべてのシグナル)、こ れが起こる可能性があります。

考えられる原因

対策

コンパクトフラッシュカードが満
 通信の問題を修正します。
 ボ
 データレートを下げます。

機器ログブック

Method		Instrument run started	09:44:46	11/20/05
1200 DAD	1	Power on	10:07:24	11/20/05
1200 DAD	1	UV-lamp on	10:07:24	11/20/05
1200 DAD	1	Vis-lamp on	10:07:24	11/20/05
1200 DAD	1	No Run data available in device!	10:07:24	11/20/05
CP Macro		Analyzing rawdata SHORT_02.D	10:07:25	11/20/05
Method		Instrument Error - Method/Sequence stopped	10:07:25	11/20/05
Method		Method aborted	10:09:52	11/20/05

図 64 機器ログブック

注記

ログブックに、通信損失(停電)は表示されません。回復(電源オン、ラン プ点灯)を示すだけです。

10 エラー情報 DAD 検出器エラー メッセージ



1220 Infinity LC

11 メンテナンス

PM の作業範囲とチェックリスト 279 注意と警告 280 溶媒送液システム 282 概要 282 溶媒フィルタのチェックとクリーニング 284 パッシブインレットバルブの交換 286 アウトレット バルブの交換 288 パージバルブフリットまたはパージ バルブの交換 290 ポンプヘッドアセンブリの取り外し 293 ポンプ シールの交換とシール馴らし作業 294 プランジャの交換 297 ポンプヘッドアセンブリの再設置 299 デュアル チャンネル グラジエントバルブ (DCGV) の交 換 300 マニュアルインジェクタ 303

メンテナンス作業の概要 303 マニュアルインジェクタのフラッシング 303 注入バルブシールの交換 304

オート サンプラ 307

概要 307

- ニードルアセンブリの交換 310 ニードルシートアセンブリの交換 315 ロータ シールの交換 317
- メタリングシールの交換 321
- グリッパアームの交換 325



11 メンテナンス

DAD 検出器エラー メッセージ

可変波長検出器 (VWD) 327 概要 327 重水素ランプの交換 328 フローセルの交換 330 フローセルの修理 332 キュベットホルダの使用 335 リークの補正 337 ダイオードアレイ検出器 (DAD) 338 メンテナンスの概要 338 モジュールのクリーニング 339 ランプの交換 340 フローセルの交換 343 フローセルのメンテナンス 346 標準フロー セルのキャピラリの交換 349 酸化ホルミウム フィルタのクリーニングまたは交 換 354 リークの補正 357 リーク処理システム部品の交換 358 HPLC システムでの藻の繁殖 359 モジュールのファームウェアの交換 361

この章では、機器のメンテナンスに関する一般情報を説明します。

メンテナンス 11

PM の作業範囲とチェックリスト

PM の作業範囲とチェックリスト

アジレントメンテナンスの作業範囲とチェックリスト

アジレントメンテナンス (PM) を実施するには、PM 作業範囲と PM チェッ クリストに段階的に従ってください。PM 作業範囲と PM チェックリストの ドキュメントは、Lab Advisor ソフトウェアの DVD に収録されています。

11 メンテナンス 注意と警告

注意と警告

警告
 電源コードが差し込まれている限り、電源を切っていても、モジュールは部分的に通電しています。
 感電やその他の身体傷害の危険性があります。モジュールの修理作業により人身障害に至る恐れがあります。たとえば、モジュールカバーが開いていて機器が電源に接続されている場合の感電などです。
 →上部カバーを取り外し、電源ケーブルを差し込んだ状態で、モジュールの調整、メンテナンス、修理を決して行わないでください。
 → 電源ケーブルコネクタの安全レバーによって、電源を接続したままモジュールのカバーを外すことはできません。カバーが取り外されている時に、電力線を決して差し込まないでください。

警告 尖った金属の先端

機器の尖った先端部分が怪我の原因になることがあります。

→ 人身障害を防ぐために、尖った金属部分に触れる際には注意して ください。

警告

有毒、可燃性および有害な溶媒、サンプル、試薬

- 溶媒、サンプル、および試薬の取り扱いには、健康や安全性を脅か す危険性が伴うことがあります。
- → これらの物質を取り扱う場合は、供給元の提供する物質の取り扱いおよび安全データシートに記載された適切な安全手順(保護眼鏡、安全手袋、および防護衣の着用など)に従ってください。
- →使用する物質の量は、分析のために必要な最小限の量に抑えてく ださい。
- → 爆発性雰囲気の中で機器を操作することはおやめください。

注意 電子ボードおよび部品は、静電気放電(ESD)に敏感です。

- ESD は電子ボードやコンポーネントを損傷を与える可能性があります。
- →ボードは端を持つようにし、電気コンポーネントには触れないようにしてください。電子ボードやコンポーネントを取り扱う時は、必ず静電気防止キット(静電気防止ストラップなど)を使用してください。

検出器光線よる目の障害

2

本製品に使用されている光学システムのランプの光を直接目で見る と、目を傷める危険があります。

→ 重水素ランプを取り外す際は、必ず光学システムのランプをオフ にしてください。

注意外部装置の安全規格

→機器に外部装置を接続する場合は、外部装置のタイプに適した安全規格 に従ってテスト、承認されたアクセサリユニットのみを使用してください。

警告

11 メンテナンス

溶媒送液システム

溶媒送液システム

概要

Agilent 1220 Infinity LC ポンプは、簡単に修理できるように設計されて います。本節で説明する手順は、ポンプをラックの所定の位置に設置した まま行えます。

プランジャシール交換やパージ バルブシール交換などの頻繁に行う修理 は、ポンプ前面から行えます。

作業	通常の実行時期	注記
溶媒フィルタのチェッ クとクリーニング	溶媒フィルタが詰まっ た場合	グラジエントパフォー マンスの問題、断続的 な圧力変動
パッシブインレットバ ルブの交換	内部でリークが発生し た場合	圧力リップルが不安定 になる場合。確認のた め、 リークテスト を実 行してください
アウトレットボールバ ルブの交換	内部でリークが発生し た場合	圧力リップルが不安定 になる場合。確認のた め、 リークテスト を実 行してください
パージバルブフリット またはパージ バルブの 交換	内部でリークが発生し た場合	バルブを閉じたとき、 廃液出口から溶媒が滴 下する。
パージバルブフリット またはパージ バルブの 交換	フリットに汚染または 詰まりが発生した場合	フリットの両側 で 10 bar を超える圧 力降下がある場合 (パージオンの状態で 5 mL/min H ₂ 0) は、フ リットが詰まっていま す。

表 44 簡単な修理手順 - 概要

メンテナンス 11

溶媒送液システム

作業	通常の実行時期	注記
ポンプ シールの交換	シールの摩耗のためポ ンプの性能が低下した 場合	ポンプヘッドの下側に リークがある。リテン ションタイムが不安定 になる。圧力リップル が不安定になる。-確認 のため、 リークテスト を実行してください。
シール馴らし作業	ポンプ シールを交換し た後	
ポンプヘッドアセンブ リの取り外し	シールまたはプラン ジャの交換前に	
プランジャの交換	傷がついた場合	シールの寿命が通常よ り短い。シールの交換 時にプランジャを点検 してください。

表 44 簡単な修理手順 - 概要

警告

警告

電源を切っていても、機器は部分的に通電しています。

フロントパネルのスイッチがオフになっている場合でも、パワー サ プライは少量の電力を使用します。

→ Agilent 1220 Infinity LC ポンプを電源から切り離すには、電源 コードのコネクタを外します。

📕 尖った金属の先端

機器の尖った先端部分が怪我の原因になることがあります。

→ 人身障害を防ぐために、尖った金属部分に触れる際には注意して ください。 11 メンテナンス 溶媒送液システム

- 警告
 キャピラリまたはチューブのフィッティングを開けると、溶媒が漏れ出す可能性があります。
 毒性や危険性のある溶媒と試薬の取り扱いには、健康上のリスクがある恐れがあります。
 >試薬メーカーから提供されている取扱説明書および安全データシートの記載に従って、適切な安全手順(ゴーグル、安全手袋、
 - 防護衣を着用するなど)を守ってください。特に、毒性の溶媒や 危険な溶媒を使用する場合は、注意してください。
- 注意 電子ボードおよび部品は、静電気放電(ESD)に敏感です。 ESD は電子ボードやコンポーネントを損傷する可能性があります。 → 損傷を避けるために、電子ボードと部品に触れる場合は、静電気防止用 保護具を必ずご使用ください。

溶媒フィルタのチェックとクリーニング

- 警告 小さな粒子がキャピラリとポンプのバルブを永久的に詰まらせる恐れがあります。 Agilent 1220 Infinity LC ポンプの損傷
 - → 溶媒は必ずろ過します。
 - → 溶媒インレットフィルタなしにポンプを決して使用しないでください。

注記

フィルタの状態が良好ならば、溶媒チューブから溶媒が自由に滴下します(静水圧)。しかし、溶媒フィルタが部分的に詰まっている場合は、溶媒チューブから溶媒はほとんど滴下しません。

メンテナンス 11

溶媒送液システム

溶媒フィルタのクリーニング

日時: 溶媒フィルタが詰まった場合

必要な部品: 説明 濃硝酸(65 %) 再蒸留水 ビーカー

必要な準備: 注入ロバルブから溶媒インレットチューブを取り外します。

- 1 詰まった溶媒フィルタをボトルヘッドアセンブリから取り外し、その フィルタを濃硝酸(65%)のビーカーに 1 時間入れておきます。
- 2 フィルタをLC クラスの水で徹底的に洗浄します(カラムの中には、濃 硝酸によって損傷するものもありますので、硝酸を完全に除去してくだ さい。pH 指示薬で確認してください)。
- 3 フィルタを元どおりに取り付けます。

11 メンテナンス 溶媒送液システム

パッシブインレットバルブの交換

- 日時: 内部でリーク(逆流)が発生した場合
- **必要なツール: 説明** レンチ、14 mm ピンセット 1 組
- **必要な部品: 部品番号 説明** G4280-60036 パッシブインレットバルブ
- **必要な準備**: ポンプの下に溶媒ボトルを置きます。
 - 1 上部前面カバーを外します。
 - インレットバルブから溶媒インレットチューブを切り離してください (静水圧により溶媒が漏れることもあるのでご注意ください)。
 - 3 インレットバルブ (オプション)からアダプタを取り外します。

溶媒送液システム

4 14 mm レンチを使用して、注入ロバルブを緩め、ポンプヘッドからバル ブを取り外します。



図 65 パッシブインレットバルブ

パッシブインレットバルブ:部品番号 G4280-60036

1 0-リング:部品番号 0905-1684

- 5 新しいバルブをポンプヘッドに差し込みます。
- 6 14 mm レンチを使用して、手で締められるだけナットを回します。
- 7 インレットバルブ (オプション) にアダプタを再び接続します。
- 8 溶媒インレットチューブをアダプタに再接続します。
- 9 前面カバーを元に戻します。

11 メンテナンス 溶媒送液システム

アウトレット バルブの交換

- 日時: 内部でリークが発生した場合
- **必要なツール: 説明** レンチ、14 mm
- 必要な部品: 部品番号 説明G1312-60067 アウトレットバルブ 1220/1260
- **必要な準備:** 主電源スイッチでポンプの電源を切ります 上部前面カバーを外します
 - 1/4 インチレンチを使用して、アウトレット バルブからバルブキャピラ リを外します。
 - 2 14 mm レンチを使用してバルブを緩め、ポンプ本体から取り外します。
 - 3 新しいバルブが正しく取り付けられたか、およびゴールドシールが付いているかを確認してください(ゴールドシールが変形している場合は、交換してください)。



- 図 66 アウトレットボールバルブ
- 4 アウトレット バルブを元どおりに取り付けて、バルブを締めます。


5 バルブキャピラリを再接続します。

パージバルブフリットまたはパージ バルブの交換

日時:フリット:プランジャシールの交換時、またはフリットが汚れているか、詰
まっている場合(フリットの両側で 10 bar を超える圧力降下が生じ、それ
はパージバルブを開いた状態で 5 mL/min の流速で H20 を送液時)

パージバルブ: 内部でリークが発生した場合

- **必要なツール: 説明** スパナ、1/4 インチ レンチ、14 mm ピンセット 1 組 または つまようじ
- 必要な部品: 番号 部品番号 説明
 1 01018-22707 PTFE フリット(5 個)
 1 G4280-60061 パージバルブ
- **必要な準備:** 主電源スイッチでポンプの電源を切ります 上部前面カバーを外します
 - 1/4インチスパナを使用して、パージ バルブからポンプのアウトレット キャピラリを切り離します。
 - 2 廃液チューブを外します。溶媒のリークに注意してください。
 - 3 14 mm レンチを使用してパージ バルブを緩め、取り外します。
 - 4 パージバルブからシールキャップを取り外します。

溶媒送液システム

5 ピンセットまたはつまようじを使用して、フリットを取り外します。



図 67 パージ バルブの部品

- 6 前図に示した向きで、新しいフリットをパージバルブに入れます。
- 7 ゴールドシール付けのキャップを元どおりに取り付けます。
- 注記 ゴールドシールが変形している場合は、元どおりに取り付ける前に交換してく ださい。
 - 8 パージバルブをポンプヘッドの中に差し込み、ポンプアウトレットキャ ピラリと廃液チューブを設置します。

溶媒送液システム



9 パージバルブを締めて、アウトレットキャピラリと廃液チューブを再接続します。

図 68 パージ バルブの交換

ポンプヘッドアセンブリの取り外し

- 日時:
 シール交換前

 プランジャ交換前
- **必要なツール: 部品番号 説明** スパナ、1/4 インチ 8710-2392 4.0 mm 六角レンチ、長さ 15 cm、T ハンドル
- **必要な準備:** 主電源スイッチでポンプの電源を切ります

警告 ポンプヘッドを取り外した状態でポンプを起動すると、ポンプドラ イブを損傷する恐れがあります。

→ポンプヘッドを取り外した状態でポンプを起動しないでください。

- 1 上部前面カバーを外します。
- 2 1/4 インチのレンチを用い、アウトレットキャピラリを取り外します。
- 3 アウトレットバルブからキャピラリを切り離します。
- 4 廃液チューブを取り外し、インレットバルブチューブを切り離します。
- 5 ポンプヘッドの下側でキャピラリを取り外します。
- 6 4 mm 六角レンチを使用して 2 個のポンプヘッドネジを徐々に緩めて取 り外し、ポンプドライブからポンプヘッドを取り外します。



11 メンテナンス 溶媒送液システム

ポンプ シールの交換とシール馴らし作業

ポンプ シールの交換

- 日時: リークテストの結果、シールのリークが発生しているとわかった場合
- 必要なツール: 部品番号 説明 スパナ、1/4 インチ
 8710-2392 4.0 mm 六角レンチ、長さ 15 cm、T ハンドル
 必要な部品: 部品番号 説明
 - 5063-6589計量シール(2個入)、100µ1分析ヘッド用0905-1420PEシール(2個入)5022-2159リストリクションキャピラリ

必要な準備: 主電源スイッチでポンプの電源を切ります。 上部フロント カバーを外します。



溶媒送液システム



シール馴らし作業

注意 次の部品ではこの手順が必要です。シール(2 個入)(5063-6589)のみ。 次の部品を損傷させます。PE シール(2 個入)(0905-1420)。

- → 順相アプリケーション用にシール馴らし作業を決して行わないでください。
- 1 100 mL のイソプロパノール入りボトルを溶媒キャビネットに設置して、 ボトルヘッドアセンブリ(チューブを含む)をボトルに取り付けます。
- 2 ボトルヘッドからのインレットチューブを直接注入ロバルブに接続します。
- 3 次の部品の一端を接続する際、リストリクションキャピラリ (5022-2159)をパージバルブに接続します。もう一方の端を廃液コンテ ナの中に挿入します。
- **4** パージバルブを開き、イソプロパノールを使って流量 2 mL/min で、シ ステムを 5 min パージします。
- 5 パージバルブを閉じて、流量を 350 bar の圧力に達するのに十分な速度 に設定します。
- 6 シールを馴らすため、この圧力で 15 min 送液します。
- ポンプをオフにし、パージ バルブをゆっくり開けて、システムから圧 力を解放します。
- 8 リストリクション キャピラリを切り離し、アプリケーション用の溶媒 を入れたボトルを元どおりに取り付けます。
- 9 システムをアプリケーションに使用する溶媒で洗浄します。

溶媒送液システム

プランジャの交換

- **日時**: プランジャに傷がついたとき
- **必要なツール: 部品番号 説明** スパナ、1/4 インチ 8710-2392 4.0 mm 六角レンチ、長さ 15 cm、T ハンドル **必要な部品: 部品番号 説明**
- 5067-4695 サファイア製ピストン
- **必要な準備**: 主電源スイッチでポンプの電源を切ります。 上部フロント カバーを外します。



溶媒送液システム



メンテナンス 11 溶媒送液システム

ポンプヘッドアセンブリの再設置

ポンプの再組み立て時 日時:

必要なツール: 説明

六角レンチ、4 mm

1 ポンプヘッドアセンブリをポンプドライブにスライドさせ、4 mm 六角 レンチを使用してポンプヘッドネジを、トルクを増しつつ、徐々に締め ます。



ポンプヘッドネジ

- 2 キャピラリとチューブを再び接続します。
- 3 前面カバーを元に戻します。

メンテナンス
 溶媒送液システム

デュアル チャンネル グラジエントバルブ (DCGV) の交換

- **必要なツール: 部品番号 説明** 8710-0899 ドライバ、Pozidriv #1
- 必要な部品: 番号 部品番号 説明
 - 1 G4280-60004 デュアルチャンネルグラジエントバルブ

必要な準備: 電源スイッチでポンプの電源を切ります。 ポンプの機構にアクセスできるように、上部フロント カバーを取り外しま す。

注記 デュアル チャンネル グラジエントバルブの寿命を維持するために、特にバッ ファ溶液を使用する場合は、定期的にバルブをフラッシングしてください。 バッファ溶液を使用する場合は、バルブのすべてのチャンネルを水でフラッ シュして、バッファ溶液から沈殿が析出するのを防止してください。塩の結晶 が発生すると、未使用のチャンネルに入り込んで塞いでしまうため、そのチャ ンネルにリークが発生することがあります。このようなリークがあると、バル ブの性能が低下します。Agilent 1220 Infinity LC ポンプにバッファ溶液と有 機溶媒を使用する際は、バッファ溶液をグラジエントバルブの下部ポートに接 続し、有機溶媒を上部ポートに接続することをお勧めします。有機チャンネル を塩溶液チャンネルの真上に設置するのがベストです(A :塩溶液、B : 有機 溶媒)。

溶媒送液システム



溶媒送液システム



マニュアルインジェクタ

マニュアルインジェクタ

メンテナンス作業の概要

表 45 メンテナンス作業の概要

作業	通常の実行時期	所要時間
インジェクタのフラッ シング	水性緩衝液または塩溶 液の使用後	5 min
インジェクションバル ブシールの交換	約 10000 ~ 20000 注入した後、またはバ ルブパフォーマンスに リークや摩耗の兆候が ある場合	10 min

マニュアルインジェクタのフラッシング

- 注意 水性緩衝液または塩溶液を使用すると、結晶を結成する可能性があります。 結晶が結成すると、インジェクションシールに傷を付ける恐れがあります。 → 水性緩衝液または塩溶液の使用後、水でバルブを必ずすすいでくださ い。
 - 1 バルブを [INJECT] ポジションに切り替えます。
 - 2 ポンプを用いて、サンプルループとシール溝を洗い流します。
 - ニードルポートクリーナー(バルブに付属)とシリンジを用いて、ニードルポートとベントキャピラリをフラッシュします。

マニュアルインジェクタ

注入バルブシールの交換

- 日時: 注入量の再現性が悪く インジェクタバルブがリークする
- **必要なツール: 説明** 六角レンチ、9/64 インチ (ツールキット内)
- **必要な部品: 部品番号 説明** 0101-1409 ローターシール、PEEK



マニュアルインジェクタ



マニュアルインジェクタ





オート サンプラ

概要

オート サンプラは、簡単に修理できるように設計されています。頻繁に行う修理は、Agilent 1220 Infinity LC 内に機器を設置したまま、機器の正面から行えます。これらの修理については、以下のセクションで説明します。

作業の概要

作業	通常の実行時期	所要時間
ニードルアセンブリの交換	ニードルが損傷や詰まりの兆候を示 す場合	15 min
シートアセンブリの交換	シートが損傷や詰まりの兆候を示す 場合	10 min
ロータ シールの交換	約 30000 ~ 40000 の注入後、ま たはバルブパフォーマンスにリーク や摩耗の兆候がある場合	30 min
メタリングシールの交換	オート サンプラの再現性から、 シールが摩耗していると判断できる 場合	30 min
グリッパアームの交換	グリッパアームに不具合がある場合	10 min

警告 正面パネルの電源スイッチを OFF にしても、電源では少量の電力が 使用されています。

> オート サンプラの修理作業により人身障害に至る恐れがあります。 たとえば、オート サンプラのカバーが開いていて、機器が電源に接 続されている場合の感電などです。

- → 電源コネクタに常にアクセスすることが可能か確認します。
- → カバーを開ける前に、機器から電源ケーブルを取り外します。
- → カバーが取り外されている間は、電源ケーブルを機器に接続しないでください。

内部の部品の交換

故障した内部の部品を交換する場合があります。これらの部品の交換には、 Agilent 1220 Infinity LC からオート サンプラを取り外す必要があり、 これらの修理は訓練を受けた要員だけが行う必要があります。

安全フラップ、フレックスボード

安全フラップとフレックスボードの交換は、アジレントの訓練を受けた サービス要員だけが行うことを強くお勧めします。

トランスポートアセンブリ部品

モータの調節やドライブベルトのテンション調整は、トランスポートアセ ンブリを正しく操作するために重要です。ドライブベルトとグリッパアセ ンブリの交換は、アジレントの訓練を受けたサービス要員だけが行うこと を強くお勧めします。トランスポートアセンブリには、他に現場で交換で きる部品はありません。その他のコンポーネント(フレックスボード、ス ピンドル、プラスチック製部品)に不具合がある場合、ユニット全体を交 換する必要があります。

オートサンプラのクリーニング

■ 感電の危険性

警告

- オートサンプラ内に液体が滴ると、感電事故を起こし、オートサン プラに損傷を及ぼす恐れがあります。
- → フィッティングを外す前には必ず、すべての溶媒ラインを排水してください。

オートサンプラのカバーは常に清潔に保つ必要があります。クリーニング は、水または中性洗剤水溶液に浸した柔らかな布を用いてクリーニングし ます。多量の水分を含んだ布はオートサンプラ中に液体を入れてしまう可 能性があるため、使用しないでください。

メンテナンス機能

特定のメンテナンス作業では、部品にアクセスしやすいように、ニードル アーム、メタリングデバイス、グリッパアセンブリを所定の位置に移動さ せる必要があります。メンテナンス機能は、これらのアセンブリを適切な メンテナンスポジションに移動させます。

11 メンテナンス オート サンプラ

ニードルアセンブリの交換

- 日時:
 ニードルが明らかに破損している場合

 ニードルが詰まっている場合
- **必要なツール: 説明** スパナ、1/4 インチ(HPLC ツールキットに付属) 2.5 mm 六角レンチ(HPLC ツールキット内) プライヤー 1 組
- **必要な部品: 番号 部品番号 説明** 1 G1313-87201 ニードルアセンブリ

必要な準備: ニードル交換(機器ユーティリティまたはLab Advisor ソフトウェアのツー ル 機能内)を選択して、開始を選択します。 ニードルがニードルシートの上、約 15 mm に位置する場合、上部フロント カバーを取り外します。

警告 人身傷害

怪我をしないように、オートサンプラの操作中はニードルエリアに 指を触れないでください。

- → 安全フラップを所定の位置から折り曲げたり、安全カバーを取り 外さないでください。
- → グリッパがニードルの下にあるときにグリッパにバイアルを挿入したり、グリッパからバイアルを取り出さないでください。





オート サンプラ



次のステップ:

9 この操作を完了するには:フロントカバーを取り付けます。

10 終了(ツール機能ニードル交換内)を選択します。

ニードルシートアセンブリの交換

 日時:
 シートが明らかに破損している場合

 シートキャピラリが詰まっている場合

必要なツール: 説明 スパナ、1/4 インチ(HPLC ツールキットに付属) マイナスドライバ

- 必要な部品:番号 部品番号説明1G1313-87101ニードルシートアセンブリ (内径 0.17 mm、2.3 µL)
- **必要な準備:**・ 開始を選択します(**ツール**機能内、LMD ソフトウェア**ニードル交換**)
 ・ 上部フロント カバーを外します。
 - ニードルアップコマンド(ニードル交換機能内)を使用して、ニードルを
 さらに 1 cm 持ち上げます。





オート サンプラ

ロータ シールの交換

- 日時: 注入量の再現性が悪く インジェクタバルブがリークする
- **必要なツール: 説明** 1/4 インチスパナ 六角レンチ、9/16 inch (ツールキット内)
- 必要な部品: 番号 部品番号 説明

 0101-1416 ローターシール (PEEK・

 必要な準備: ・ 上部前面カバーを取り外します。

 リークチューブを取り外します (必要な場合)。

 注意 ステータヘッドの取り外し

 ステータへッドの取り外し
 ステータフェースは、ステータへッドによって固定されています。ステータヘッドを取り外す場合は、ステータフェースがバルブの外に落ちる恐れがあります。
 - → ステータフェースへの損傷を防ぐように、バルブを慎重に取り扱います。



オート サンプラ



次のステップ:

9 リークトレイ内の廃液ホルダに廃液チューブを差し込みます。

10この操作を完了するには:フロントカバーを取り付けます。

オート サンプラ

メタリングシールの交換

- 日時: 注入量の再現性が悪く メタリングデバイスのリーク
- 必要なツール: 部品番号 説明
 スパナ、1/4 インチ (HPLC ツールキットに付属)
 4 mm 六角レンチ (HPLC ツールキット内)
 8710-2411 3 mm 六角レンチ (HPLC ツールキット内)
- 必要な部品:番号 部品番号説明15063-6589計量シール(2個入)、100µ1分析ヘッド用15063-6586ピストン

必要な準備:・ 開始(ツール機能内、LMD ソフトウェアピストン交換)を選択します。 ・ 上部フロント カバーを外します。





オート サンプラ




オート サンプラ

グリッパアームの交換

- 日時: グリッパアームの故障
- **必要なツール: 説明** まっすぐにした紙クリップ
- 必要な部品:番号 部品番号説明1G1313-60010グリッパアセンブリ
- **必要な準備:**・開始(ツール機能内、LMD ソフトウェアグリッパの交換)を選択します。
 - ・ 機器への電源を切ります。
 - 上部フロント カバーを外します。



11 メンテナンス オート サンプラ



可変波長検出器(VWD)

可変波長検出器 (WWD)

概要

このセクションでは、メインカバーを開けずに行える検出器の簡単なメン テナンスや修理手順を説明します。

表 46 検出器のメンテナンスと修理

作業	通常の実行時期	注記
重水素ランプの交換	ノイズやドリフトがアプリケーション のリミット値を超えた場合、またはラ ンプが点火しない場合。	交換後に VWD テストを行う 必要がある。
フローセルの交換	アプリケーションが、異なるタイプの フローセルを必要とする場合。	交換後に VWD テストを行う 必要がある。
フローセルの修理	リークがある、またはフローセル窓の 汚れのために強度が低下した場合。	修理後、圧力気密性テスト を実施。
リーク センサの乾燥	リークが発生した場合	リークをチェックします。

可変波長検出器(VWD)

重水素ランプの交換

- **日時**: ノイズまたはドリフトが使用目的のリミット値を超えている場合、またはランプが点灯しない場合
- **必要なツール: 説明** ドライバ、Pozidriv #1 PT3
- 必要な部品:番号部品番号説明1G1314-60100重水素ランプ
- **必要な準備**: ランプを OFF にします。

警告 高温のランプを触れることによる怪我 検出器を使用していた場合は、ランプが熱くなっています。 → その場合は、ランプが冷えるまで待ちます。



可変波長検出器(VWD)



可変波長検出器(VWD)

フローセルの交換

日時: アプリケーションが、異なるタイプのフローセルを必要とする場合、または フローセルのメンテナンスが必要な場合

必要なツール: 説明 レンチ、1/4 inch キャピラリ接続用

必要な準備: ランプを OFF にします。

1 リリース ボタンを押して下部フロント カ	2 インレットキャピラリとアウトレットキャ
バーなんし、フローセル領域にアクセフト	
ハーを外し、ノロービル領域にノクビスし	レノリゼハしより。
ます。	
	A start
	~

可変波長検出器(VWD)



可変波長検出器 (VWD)

フローセルの修理

必要な部品: 部品番号 説明
 G1314-60086 標準フローセル、10 mm、14 µL、40 bar
 G1314-65061 セルリペアキット、内容:2 x ガスケット #1、2 x ガスケット
 ト #2、2 x ウィンドウクォーツ

→ ウィンドウを取り外すためにピンセットは使用しないでください。



図 69 標準フロー セル 10 mm / 14 µL

1 リリース ボタンを押して下部フロント カバーを外し、フローセル領域 にアクセスします。

可変波長検出器(VWD)

- **2** インレットキャピラリとアウトレットキャピラリを外します。
- **3** つまみネジを同時に緩めてフローセルを外します。
- 4 フローセルの分解
 - a 4 mm 六角レンチを使用して、セルネジを外します。
 - **b** ピンセットを使用して、SST リングを外します。
 - c 粘着テープを使用して、PEEK リング、ウィンドウ、ガスケットを取 り外します。
 - d もう一方のウィンドウについても『ステップ 1 ページ 332』スルー 『ステップ 3 ページ 333』を繰り返します。部品は分けて置きます。 そうしないと混ざってしまいます。
- 5 フローセル部品のクリーニング
 - a イソプロパノールをセル穴に注ぎ、糸くずが出ない布できれいに拭き ます。
 - b エタノールかメタノールでウィンドウをきれいにします。糸くずが出ない布でウィンドウを乾かします。
- 6 フローセル再組み立て
 - a フローセルのカセットを水平に保ち、ガスケットを正しい位置に置き ます。ガスケットの穴を通して両方のセル穴が見えることを確認しま す。
- 必ず新しいガスケットを使用してください。
 - **b** ウィンドウをガスケットの上に置きます。
 - **c** PEEK リングをウィンドウの上に置きます。
 - d コニカルスプリングを挿入します。円錐スプリングがウィンドウの方 に向いていることを確認してください。逆の場合、セルネジを締める と、ウィンドウが壊れることがあります。
 - e セルネジをフローセルの中に入れて、ネジを締めます。
 - 7 セルのもう一方の側にこの作業を繰り返します。
 - 8 インレットキャピラリとアウトレットキャピラリを再接続します。

注記

可変波長検出器 (VWD)

- 9 フローセルにリークがないかテストします。リークがない場合は、フ ローセルを検出器に挿入します。
- 10 波長キャリブレーションを行い、フローセルの位置が正しいか確認しま す(『「波長ベリフィケーション/キャリブレーション」177 ページ』)。
- 11 下部フロント カバーを取り付けます。

可変波長検出器(VWD)

キュベットホルダの使用

このキュベットホルダを、フローセルの代わりに UV-Vis 検出器に取り付けることができます。スタンダードを入れた標準キュベット(例えば、米国連邦標準技術局(NIST)ホルミウムオキサイドスタンダード溶液)をその中に固定することができます。

これは、波長検証に使用することができます。

- 日時: 独自の標準を使用して装置をチェックアウトする場合
- **必要なツール: 説明** なし
- **必要な部品: 部品番号 説明** G1314-60200 キュベットホルダ 「標準サンプル」(NIST 承認酸化ホルミウムサンプルなど)入 りのキュベット

必要な準備: フローセルを外します。 標準を入れたキュベットを使用できるようにします。



可変波長検出器(VWD)



可変波長検出器(VWD)

リークの補正

- **日時**: フローセル領域またはキャピラリの接続部にリークが発生した場合
- **必要なツール: 説明** ティッシュペーパー レンチ、1/4 inch キャピラリ接続用

必要な部品: 説明

なし

- 1 下部前面カバーを外します。
- 2 ティッシュペーパを使用して、リークセンサ領域を拭いて乾かします。
- **3** キャピラリ接続部とフローセル領域にリークがないか確認し、必要な場合は処置を施します。
- 4 フロント カバーを取り付けます。



ダイオードアレイ検出器 (DAD)

ダイオードアレイ検出器 (DAD)

メンテナンスの概要

以下のページでは、メインカバーを開けずに行える検出器のメンテナンス (簡単な修理)を説明します。

表 47 メンテナンスの概要

作業	通常の実行時期	注記
モジュールの クリーニング	必要な場合	
重水素ランプまた はタングステン ランプの交換	ノイズやドリフトがアプリケーションのリ ミット値を超えた場合、またはランプが点 火しない場合。	交換後に VWD テストを行う必 要があります。
フローセルの交換	アプリケーションが、異なるタイプのフ ローセルを必要とする場合。	交換後、ホルミウムテストま たは波長キャリブレーション テストを行います。
フローセルの部品 のクリーニング または交換	リークがある、またはフローセル窓の汚れ のために強度が低下した場合。	修理後、圧力気密性テストを 実施。
酸化ホルミウム フィルタのクリー ニングまたは交換	汚れた場合。	交換後、ホルミウムテストま たは波長キャリブレーション テストを行います。
リーク センサの 乾燥	リークが発生した場合	リークをチェックします。
リーク処理 システムの交換	破損または腐蝕した場合	リークをチェックします。

ダイオードアレイ検出器(DAD)

モジュールのクリーニング

モジュールのケースは、清潔に保つ必要があります。クリーニングする際 は、少量の水または弱い洗剤を水で薄めた溶液に浸した柔らかい布を使用 してください。大量の液体を含んだ布を使用すると、液体がモジュール内 に滴下することがあります。

→ クリーニング中は多量の水分を含んだ布を使用しないでください。

→ 流路内の連結部を開く前には必ず、すべての溶媒ラインを排水してください。

11 メンテナンス ダイオードアレイ検出器 (DAD)

ランプの交換

警告

日時: ノイズまたはドリフトがアプリケーションのリミット値を超えている場合、 またはランプが点火しない場合

必要なツール: 説明 ドライバ、Pozidriv #1 PT3

- **必要な部品: 番号 部品番号 説明** 1 2140-0820 長寿命重水素ランプ「C」(黒色カバーおよび ID タグ付き)
- または 1 G1103-60001 タングステンランプ

学 おりには 男される

本製品に使用されている重水素ランプの光を直接目で見ると、 目をいためる危険があります。

→ 必ず重水素ランプを OFF にしてから、重水素ランプを外す作業を 行ってください。

警告
 高温のランプを触れることによる怪我
 検出器を使用していた場合は、ランプが熱くなっています。
 → その場合は、ランプが冷えるまで待ちます。

ダイオードアレイ検出器(DAD)



ダイオードアレイ検出器(DAD)

次のステップ:

- 5 フロント カバーを閉めます。
- 6 ユーザーインタフェースのドキュメントに記載されているように、ランプカウンタをリセットします (I.D. タグ付きのランプはリセットできません)。
- **7** ランプをオンにして、ランプを10分ウォームアップします。
- 8 『「波長ベリフィケーションおよびリキャリブレーション」199 ページ』または 『「ホルミウム オキサイドテスト」190 ページ』を実行して、UW ランプの位置が正しいか確認します。
- 9 『「強度テスト」187 ページ』を実行します。

ダイオードアレイ検出器(DAD)

フローセルの交換

日時: アプリケーションが、異なるタイプのフローセルを必要とする場合、または フローセルの修理が必要な場合

必要なツール: 説明

レンチ、1/4 inch キャピラリ接続用

必要な部品:	番号	部品番号	説明
	1	G1315-60022	標準フローセル、10 mm、13 μL、120 bar (12 MPa)
または	1	G1315-60025	セミマイクロフローセル、6 mm、5 μL、120 bar (12 MPa)
または	1	G1315-60024	マイクロフローセル、3 mm、2 μL、120 bar (12 MPa)
または	1	G1315-60015	耐高圧フローセル、6 mm、1.7 μL、400 bar (40 MPa)
または	1	G1315-68716	ナノフローセルキット、10 mm、80 nL、5 MPa

必要な準備: ランプを消灯します。

フロント カバーを取り外します。

1 フロント カバーを開き、フローセル領域に	2 フローセルカバーを開きます。
アクセスします。	注記 システム設定によって、インレットキャピ ラリがモジュールから直接セルの上または 下に送られ、キャピラリ ホルダに送られな

ダイオードアレイ検出器 (DAD)



メンテナンス 11 ダイオードアレイ検出器(DAD)

7 フローセルキャピラリをユニオン ホルダに	次のステップ:
挿入します(上部が注入口、下部が出口)。	
つまみネジを締めて、廃液チューブ(下部)	8 『「波長ベリフィケーションおよびリキャリ
をユニオンに再接続します。	ブレーション」199 ページ』または 『「ホル
注記 リークをチェックするには、溶媒を流して フローセル (セルコンパートメントの外側) とすべてのキャピラリ接続部を確認します。	ミウムオキサイドテスト」190 ページ』を実 行し、フローセルの位置が正しいか確認し ます。 9 フロント カバーを閉じます。
注入口 出口 つまみネジ 廃液	

ダイオードアレイ検出器 (DAD)

フローセルのメンテナンス

日時: フローセルにリークまたは汚れ(光スループットが低下している)があり、 修理が必要な場合

必要なツール: 説明 レンチ、1/4 inch キャピラリ接続用 六角レンチ、4 mm つまようじ

必要な部品: 説明 部品については、『「標準フロー セル」388 ページ』を参照してください。

必要な準備: 流量をオフにします。 フロント カバーを取り外します。 フローセルを外します(『「フローセルの交換」343 ページ』を参照)。

ダイオードアレイ検出器(DAD)



ダイオードアレイ検出器(DAD)

次のステップ:

- **3** 4 mm 六角レンチを使用して、ウィンドウネジを締め付けます(手でいっぱいに締め付けた後、 レンチを使用してさらに 4 分の 1 回転締める)。
- 4 キャピラリを再接続します。『「フローセルの交換」343ページ』を参照。
- 5 リークテストを実行します。
- 6 フローセルを挿入します。
- 7 前面カバーを取り付けます。
- 8 『「波長ベリフィケーションおよびリキャリブレーション」199 ページ』または 『「ホルミウム オキサイドテスト」190 ページ』を実行し、フローセルの位置が正しいか確認します。

ダイオードアレイ検出器(DAD)

標準フロー セルのキャピラリの交換

日時: キャピラリがつまった場合

必要なツール: 説明 レンチ、1/4 inch キャピラリ接続用 レンチ、4 mm (キャピラリ接続用) ドライバ、Pozidriv #1 PT3

必要な部品: 説明 部品については、『「標準フロー セル」388ページ』を参照してください

必要な準備: ランプを消灯します。 フロント カバーを取り外します。 フローセルを外します(『「フローセルの交換」343 ページ』を参照)。

注記 この作業の説明はすべて、セルのデフォルト オリエンテーション(製造時) に基づいています。両方のキャピラリを下部または上部(キャピラリからカラ ムまでの経路に応じて)に送られるように、配管熱交換器/キャピラリとセル ボディを鏡面対称に固定できます。

注記 フローセルボディのフィッティングは、低デッド ボリューム用の特別なタイ プで、他のフィッティングに適合しません。 フィッティングを再度締める場合は、注意して締めるようにします(手できっ ちり締め、さらにレンチで 1/4 回転)。さもなければ、フローセルが損傷また は詰まることがあります。

ダイオードアレイ検出器 (DAD)



ダイオードアレイ検出器(DAD)



ダイオードアレイ検出器 (DAD)



メンテナンス 11 ダイオードアレイ検出器 (DAD)

13 熱交換器をクランプ ユニットに固定し、フ	14 インレットキャピラリをフローセルボディ
ローセルホティを熱父換畚に固正します。	に固定するにのに、ノイツテインクをよう 手でいっぱいに締め付けます。次に、4 mm
	レンチを使用してさらに 4 分の 1 回転締め
	ます。
	O A
A. Office	
15 穴に対してホルダが中心にあるか確認しま す。必要に応じて、ホルダのネジで調整し	次のステップ:
15 穴に対してホルダが中心にあるか確認します。必要に応じて、ホルダのネジで調整します。	次のステップ: 16 キャピラリを再接続します。『「フローセル
15 穴に対してホルダが中心にあるか確認します。必要に応じて、ホルダのネジで調整します。	次のステップ: 16 キャピラリを再接続します。『「フローセル の交換」343 ページ』を参照。
15 穴に対してホルダが中心にあるか確認します。必要に応じて、ホルダのネジで調整します。	 次のステップ: 16 キャピラリを再接続します。『「フローセルの交換」343 ページ』を参照。 17 リークテストを実行します。
15 穴に対してホルダが中心にあるか確認しま す。必要に応じて、ホルダのネジで調整し ます。	 次のステップ: 16 キャピラリを再接続します。『「フローセルの交換」343 ページ』を参照。 17 リークテストを実行します。 18 フローセルを挿入します。
15 穴に対してホルダが中心にあるか確認します。必要に応じて、ホルダのネジで調整します。	 次のステップ: 16 キャピラリを再接続します。『「フローセルの交換」343ページ』を参照。 17 リークテストを実行します。 18 フローセルを挿入します。 19 前面カバーを取り付けます。
15 穴に対してホルダが中心にあるか確認しま す。必要に応じて、ホルダのネジで調整し ます。	 次のステップ: 16 キャピラリを再接続します。『「フローセルの交換」343ページ』を参照。 17 リークテストを実行します。 18 フローセルを挿入します。 19 前面カバーを取り付けます。 20 『「波長ベリフィケーションおよびリキャリ
15 穴に対してホルダが中心にあるか確認します。必要に応じて、ホルダのネジで調整します。	 次のステップ: 16 キャピラリを再接続します。『「フローセルの交換」343ページ』を参照。 17 リークテストを実行します。 18 フローセルを挿入します。 19 前面カバーを取り付けます。 20 『「波長ベリフィケーションおよびリキャリブレーション」199ページ』または『「ホルミウムオキサイドテスト」190ページ』を実
15 穴に対してホルダが中心にあるか確認しま す。必要に応じて、ホルダのネジで調整し ます。	 次のステップ: 16 キャピラリを再接続します。『「フローセルの交換」343ページ』を参照。 17 リークテストを実行します。 18 フローセルを挿入します。 19 前面カバーを取り付けます。 20 『「波長ベリフィケーションおよびリキャリブレーション」199ページ』または『「ホルミウムオキサイドテスト」190ページ』を実行し、フローセルの位置が正しいか確認し
15 穴に対してホルダが中心にあるか確認しま す。必要に応じて、ホルダのネジで調整し ます。	 次のステップ: 16 キャピラリを再接続します。『「フローセルの交換」343ページ』を参照。 17 リークテストを実行します。 18 フローセルを挿入します。 19 前面カバーを取り付けます。 20 『「波長ベリフィケーションおよびリキャリブレーション」199ページ』または『「ホルミウムオキサイドテスト」190ページ』を実行し、フローセルの位置が正しいか確認します。
15 穴に対してホルダが中心にあるか確認します。必要に応じて、ホルダのネジで調整します。	 次のステップ: 16 キャピラリを再接続します。『「フローセルの交換」343ページ』を参照。 17 リークテストを実行します。 18 フローセルを挿入します。 19 前面カバーを取り付けます。 20 『「波長ベリフィケーションおよびリキャリブレーション」199ページ』または『「ホルミウムオキサイドテスト」190ページ』を実行し、フローセルの位置が正しいか確認します。

1220 Infinity LC

ダイオードアレイ検出器(DAD)

酸化ホルミウム フィルタのクリーニングまたは交換

日時: ホルミウムオキサイドフィルタが汚れた場合

必要なツール: 説明 ドライバ、Pozidriv #1 PT3 マイナスドライバー レンチ、1/4 inch キャピラリ接続用 ピンセット 1 組

必要な部品: 番号 部品番号 説明 1 79880-22711 ホルミウムオキサイドフィルター

必要な準備: ランプを消灯します。 フロント カバーを取り外します。 フローセルを外します(『「フローセルの交換」343 ページ』を参照)。

注記 『「ホルミウムフィルタの波長の証明書」417ページ』も参照してください。

ガラスは、標準の環境条件下でも表面に膜を形成する傾向があります、これは 他のいくつかのガラス表面でも見られ可能性の有る現象で、ガラスの組成と何 らかの関係があります。膜が装置に影響を及ぼすという指摘はありません。光 を著しく散乱する薄膜の場合でも、ピーク位置のシフトは予想されません。吸 光度にわずかな変化が起こるかもしれません。光路中の他のコンポーネント (レンズ、ウィンドウなど)も時間とともにその動きを変えていきます。

ダイオードアレイ検出器 (DAD)



ダイオードアレイ検出器 (DAD)



ダイオードアレイ検出器(DAD)

リークの補正

- **日時**: フローセル領域、熱交換器、またはキャピラリ接続部でリークが発生した場合
- **必要なツール: 説明** ティッシュペーパー レンチ、1/4 inch キャピラリ接続用
- **必要な準備:** 前面カバーを外します。
 - ティッシュペーパを使用して、リークセンサ箇所とリーク受けを拭いて 乾かします。
 - 2 キャピラリ接続部とフローセル領域にリークがないか確認し、必要な場合は補正します。



図 70 リークがないか確認します。

3 前面カバーを取り付けます。

ダイオードアレイ検出器 (DAD)

リーク処理システム部品の交換

- 日時: 部品が腐食したか破損した場合
- **必要なツール**: なし

必要な部品:

- 番号 部品番号 説明
 - 1 5061-3388 漏斗
 - 1 5041-8389 漏斗ホルダ
 - 5062-2463 フレックス チューブ、ポリプロピレン製、内径
 6.5 mm、5 m
- **必要な準備:** 前面カバーを外します。
 - 1 漏斗を漏斗ホルダから外します。
 - 2 漏斗をチューブとともに外します。
 - 3 漏斗をチューブとともに正しい位置に挿入します。
 - 4 漏斗を漏斗ホルダに挿入します。



図 71 リーク処理システム部品の交換

5 前面カバーを取り付けます。

HPLC システムでの藻の繁殖

HPLC システムでの藻の繁殖

HPLC システム内に藻が存在するとさまざまな問題が生じ、誤って機器やア プリケーションの問題として診断されることがあります。藻は水系溶媒中 で、特に 4 - 8 の pH 範囲で繁殖します。藻の繁殖は燐酸や酢酸など、 緩衝塩の存在で加速されます。藻は光合成によって成長するため、光もま たその繁殖を促進します。蒸留水であっても、しばらくすると小さな藻が 繁殖します。

藻が関連した機器の問題

藻は、HPLC システムのどこにでも沈着、成長し、次の問題を引き起こしま す。

- ボールバルブ、注入口あるいは出口に沈着し、流量が不安定となるか、 ポンプ全体が故障します。
- 孔径の小さな溶媒インレットフィルタを塞ぎ、流量が不安定となるか、 またはポンプ全体が故障します。
- ・ 通常インジェクタの前にある小さな孔径の高圧溶媒フィルタを塞ぎ、シ ステム圧力が上昇します。
- カラムの前に付いているフィルタを塞ぎ、システム圧力が上昇します。
- 検出器のフローセル窓が汚れ、ノイズレベルが高くなります。検出器は 流路内の最後のモジュールであることから、この問題はあまり見られま せん。

HPLC システムでの藻の繁殖

Agilent 1220 Infinity LC で見られる現象

ヘリウム脱気を利用した HP 1090 や HP 1050 シリーズなどの他の HPLC シ ステムとは対照的に、脱気にヘリウムを用いていない、Agilent 1220 Infinity LC などのシステムでは、藻が繁殖する可能性が高くなります (ほとんどの藻は繁殖に酸素と光が必要です)。

Agilent 1220 Infinity LC に藻が存在すると、次のような問題が生じる可 能性があります。

- 次の詰まり PTFE フリット(5 個)(01018-22707)(パージバルブアセンブリ)およびカラムフィルタ、システム圧力が上昇します。藻は、フィルタ上に白色または黄色がかった白色の沈着物として現れます。通常、ピストンシールの標準的な摩耗による黒い粒子は、PTFE フィルタを短期間の使用で詰まらせることはありません。
- 溶媒フィルタ(ボトルヘッドアセンブリ)の寿命が短くなる。溶媒ボトル内の溶媒フィルタが詰まった場合、特に部分的に閉塞した場合は原因の発見が困難で、グラジエント性能の低下や断続的な圧力変動として現れることがあります。
- ・藻の繁殖は、ボールバルブなど、溶媒の流路にあるコンポーネントの故
 にの原因となることもあります。

藻の発生の予防 / 軽減

- 必ず新しく調整した溶媒、特に、約 0.2 µm のフィルタでろ過した脱塩 水を用いてください。
- 移動相を流さずに、数日間、装置に放置することは避けてください。
- 古い移動相は必ず廃棄してください。
- ・以下を使用します。溶媒ボトル、褐色(9301-1450) 水性移動相用、装置に標準付属 -
- 可能であれば、数 mg/L のアジ化ナトリウム、あるいは数パーセントの 有機溶媒を水性移動相に加えてください。
メンテナンス 11

モジュールのファームウェアの交換

モジュールのファームウェアの交換

新しいファームウェアをインストールする必要がある場合

- 新しいバージョンにより、古いバージョンの問題を解決する場合
- すべてのシステムを同じ(バリデーション済み)リビジョンに保つため

古いファームウェアをインストールする必要がある場合

- すべてのシステムを同じ(バリデーション済み)リビジョンに保つため、
 または
- 新しいファームウェアの新しいモジュールをシステムに追加する場合
- サードパーティ製ソフトウェア用に特別なバージョンが必要な場合。

必要なツール: 説明

日時:

LAN/RS-232 ファームウェア更新ツール

- または Agilent 診断用ソフトウェア
- または インスタントパイロット G4208A (モジュールがサポートしている場合のみ)

必要な部品: 番号 説明

1 Agilent ホームページからのファームウェア、ツール、およびドキュメ ント

必要な準備: ファームウェア更新ツールに付属するドキュメントをお読みください。

モジュールのファームウェアをアップグレード / ダウングレードするに は、次の操作を行います。

- 必要なモジュールファームウェア、最新の LAN/RS-232 ファームウェア 更新ツール、アジレントウェブサイトにある付属文書をダウンロードし ます。
 - http://www.chem.agilent.com/scripts/cag_firmware.asp.
- 2 モジュールにファームウェアを読み込むには、付属のドキュメントの手順に従います。

モジュール特定情報

このモジュールの特定情報はありません。

1220 Infinity LC

11 メンテナンス モジュールのファームウェアの交換



1220 Infinity LC

12 メンテナンス用部品

1220 Infinity LC システム 364 システム部品 364 ヒューズ 365 溶媒送液システム 366 注入システム 375 マニュアルインジェクタ 375 オート サンプラ 377 カラムオーブン 383 検出器 384 可変波長検出器(VWD) 384 ダイオードアレイ検出器(DAD) 388

この章では、メンテナンス用部品について説明します。



12 メンテナンス用部品

1220 Infinity LC システム

1220 Infinity LC システム

システム部品

システム部品番号

部品番号	説明
0950-4997	電源
G4280-65050	Agilent 1220 Infinity LC メインボード
G4280-68713	キャビネットキット
G4280-60102	フロントドア、上部
G4280-60001	フロントドア、下部
G4280-65001	電源スイッチボード
G4280-65802	FSL ボード(ステータス LED ボード)
5067-5378	接続チューブ、DCGV から PIV
G4280-68708	SSV
G4280-80004	ファン
8121-1833	電源スイッチケーブル
G4280-81602	ケーブル、ステータス LED
G4280-81620	温度センサーケーブル
G4280-40007	ライトパイプ、ステータス
G4280-40016	ライトパイプ、ショート
G4280-44013	リークプレーン、マニュアルインジェクション
G4280-44500	リークパネル、下部
G4280-44501	リークプレーン、ポンプ用
G4280-44502	リークプレーン、ALS

メンテナンス用部品 12

1220 Infinity LC システム

部品番号	説明
G4280-44016	ホルダ、温度センサー
5061-3356	リークセンサー

ヒューズ

5 メインボードヒューズ

部品番号	説明
2110-1417	ヒューズ 3.15 A250 V

注記 各ヒューズの隣には LED があります。赤色 LED はヒューズが切れていること を示します。ヒューズの 1 つが切れると、電源スイッチの緑 LED が点滅しま す。

- ヒューズ F1 (デガッサ、ポンプ、インジェクタモータ)
- ・ ヒューズ F2 (インジェクタセンサ、カラムオーブン、拡張 24 V コネクタ)
- ヒューズ F3 (プロセッサコア、+5 V、+15 V、-15 V のメインボードで 供給)
- ヒューズ F4 (VWD、D2- ランプを含む)
- ヒューズ F5 (VWD ヒーター、ファン)
- 1 ヒューズ ネットフィルタ

部品番号 説明

2110-1004 ヒューズ 10 A t (2x)

12 メンテナンス用部品 溶媒送液システム

浴妹这次システム

溶媒送液システム

ポンプヘッドアセンブリ(シール ウォッシュなし)

品目		部品番号	説明
		G1312-60056	シールウォッシュのないポンプヘッド 1200 SL
	1	5063-6586	サファイア製ピストン
	2	G1311-60002	ピストンハウジング
	3	5067-1560	サポートリング SL(シールウォッシュなし)
	4	5062-2484	ガスケット、ウォッシュ用シール用(6 個入)
	5	5042-8952	シールホルダ
	6	5063-6589	ピストンシール PTFE、カーボン混合、黒(2 個 入)、デフォルト部品
または		0905-1420	PE シール(2 個入)
	7	G1311-25200	ポンプチャンバハウジング
	8	G1312-60066	パッシブインレットバルブ 1220/1260
	9	G1312-60067	アウトレットバルブ 1220/1260
	10	5042-1303	ロックネジ
	11	G4280-60061	パージバルブ
	12	0515-2118	ポンプヘッドネジ(M5、60 mm)

品目 シールウォッシュのないポンプヘッド 1200 SL (G1312-60056) 1 ~ 7、10、12 を含む

メンテナンス用部品 12 溶媒送液システム



図 72 ポンプヘッドアセンブリ(シールウォッシュオプションなし)

ポンプヘッドアセンブリ (シール ウォッシュ オプション付き)

品目		部品番号	説明
		G1312-60045	ポンプヘッドアセンブリ(シールウォッシュ 付き)
	1	5063-6586	サファイア製ピストン
	2	G1311-60002	ピストンハウジング
	3	01018-60027	サポートリングのシールウォッシュ
	4	0905-1175	ウォッシュシール (PTFE)
または		0905-1718	ウォッシュ用シール PE
		0890-1764	チューブ(シールウォッシュ)
	5	5062-2484	ガスケット、ウォッシュ用シール用(6 個入)
	6	5042-8952	シールホルダ
	7	5063-6589	ピストンシール PTFE、カーボン混合、黒(2 個入)、デフォルト部品
		0905-1420	PE シール(2 個入)
	8	G1311-25200	ポンプチャンバハウジング
	9	G1312-60066	パッシブインレットバルブ 1220/1260
	10	G1312-60067	アウトレットバルブ 1220/1260
	11	5042-1303	ロックネジ
	12	G4280-60061	パージバルブ
	13	0515-2118	ポンプヘッドネジ (M5、60 mm)
		01018-23702	ポンプシール取り外し用工具

品目 ポンプヘッドアセンブリ(シールウォッシュ付き)(G1312-60045)1 ~ 8、11、13 を含む アクティブ シール ウォッシュは、1220 Infinity LC 用にサポートされていま せん。連続洗浄のみです。



図 73 ポンプヘッド (シール ウォッシュ オプション付き)

注記

アウトレットボールバルブ アセンブリ

部品番号

説明

G1312-60067 アウトレットバルブ 1220/1260



パージバルブアセンブリ

品目	部品番号	説明
	G4280-60061	パージバルブ
1		バルブ本体
2	01018-22707	PTFE フリット(5 個)
3	5067-4728	シールキャップアセンブリ



パッシブインレットバルブ アセンブリ

部品番号

説明

G1312-60066 パッシブインレットバルブ 1220/1260



ボトル ヘッド アセンブリ

品目	部品番号	説明
	G1311-60003	ボトルヘッドアセンブリ
1	5063-6598	ロックリング付きフェラル(10 個入)
2	5063-6599	チューブネジ(10 本入)
3		ワイヤマーカ
4	5062-2483	溶媒チューブ、5 m
5	5062-8517	フリットアダプタ(4 個入)
6	5041-2168	溶媒インレットフィルタ、孔径 20 μm



12 メンテナンス用部品 溶媒送液システム

配管

品目	部品番号	説明
1	G1312-67305	アウトレットキャピラリ、ポンプ - 注入装置
	G1311-60003	ボトルヘッドアセンブリ、ボトルからパッシブイ ンレットバルブ またはデガッサまで
2	G4280-60034	溶媒チューブ、デガッサから DCGV まで
3	G4280-81300	キャピラリ、第 1 プランジャからダンパーまで
4	G4280-81301	キャピラリ、ダンパーから第 2 プランジャまで
5	5067-5378	接続チューブ、DCGV から PIV
6	5062-2461	廃液チューブ、5 m (再注文パック)
	G1311-60065	ダンパアイソクラティック / クォータナリポンプ 600 bar
	G1311-60001	ポンプドライブ
	G4280-60004	デュアルチャンネルグラジエントバルブ
	3160-1017	ファン



注入システム

マニュアルインジェクタ

インジェクションバルブアセンブリ

品目	部品番号	説明
	5067-4202	マニュアルインジェクションバルブ 600 bar、 ループとニードルポートを含む一式
	1535-4045	アイソレーションシール
2	0101-1409	ローターシール、PEEK
3	0101-1417	ステータヘッド
4	5067-1581	ニードルポート
5	5068-0018	ステータネジ
	8710-0060	六角レンチ、9/64 インチ



12 メンテナンス用部品 注入システム

サンプルループ

サンプル ループ ステンレス スチール

部品番号	説明
0101-1248	サンプルループ 5 μL
0100-1923	サンプルループ 10 μL
0100-1922	サンプルループ 20 µL
0100-1924	サンプルループ 50 μL
0100-1921	サンプルループ 100 µL
0101-1247	サンプルループ 200 µL
0101-1246	サンプルループ 500 µL
0101-1245	サンプルループ 1 mL
0101-1244	サンプルループ 2 mL

サンプル ループ PEEK

部品番号 説明 サンプルループ 5 μ L 0101-1241 サンプルループ 10 µL 0101-1240 0101-1239 サンプルループ 20 µL サンプルループ 50 µL 0101-1238 サンプルループ 100 µL 0101-1242 サンプルループ 200 µL 0101-1227 0101-1236 サンプルループ 500 µL サンプルループ 1 mL 0101-1235 サンプルループ 2 mL 0101-1234

オート サンプラ

オート サンプラのメインアセンブリ

部品番号	説明
G4280-60230	オートサンプラー式
G1329-60009	トランスポートアセンブリ
G4280-60027	サンプリングユニットアセンブリ(インジェクション バルブと分析ヘッドは除く)
01078-60003	アナリティカルヘッドアセンブリ、100 μL
0101-1422	インジェクションバルブ
G1313-44510	バイアルトレイ
G1313-60010	グリッパアセンブリ
G4280-87304	廃液キャピラリ
G4280-81615	ケーブル、サンプリングユニット
G4280-81616	ケーブル、サンプルトランスポート
5067-1581	ニードルポート

12 メンテナンス用部品 注入システム

アナリティカルヘッド アセンブリ

品目	部品番号	説明
	01078-60003	アナリティカルヘッドアセンブリ、100 μL
1	5063-6586	ピストン
2	0515-0850	ネジ、M4、長さ 40 mm
3	01078-23202	アダプタ
4	5001-3739	サポートシールアセンブリ
5	5063-6589	計量シール (2 個入)、100 µ1 分析ヘッド用
6	01078-27710	ヘッド本体
	0515-2118	ネジ M 5、長さ 60 mm、アセンブリ取り付け用

メンテナンス用部品 12 注入システム



12 メンテナンス用部品 注入システム

サンプリング ユニット アセンブリ

品目	部品番号	説明
	G4280-60027	サンプリングユニットアセンブリ(インジェク ションバルブと分析ヘッドは除く)
1	G1313-66503	SUD ボード
2	1500-0697	ベルトギア、計量ユニットとニードルアーム用
3	5062-8590	ステッパモータ、メタリングユニットとニードル アーム用
4	01078-87302	ループキャピラリ(100 μL)
5	01078-60003	アナリティカルヘッドアセンブリ、100 μL
6	G1313-87301	キャピラリ、インジェクションバルブから分析 ヘッドまで(160 mm × 0.25 mm)
7	G1329-44115	安全カバー
8	0101-1422	インジェクションバルブ
9	G1313-87300	廃液チューブインジェクションバルブアセンブリ (120 mm)
11	G1313-87101	ニードルシートアセンブリ (内径 0.17 mm、 2.3 µ L)
12	G1313-43204	シートアダプタ
13	G1313-44106	安全フラップ
14	G1313-68715	フレックスボード
15	G1313-87201	ニードルアセンブリ
	G1313-68713	クランプキット(ニードルクランプと 2 x クラ ンプネジを含む)

メンテナンス用部品 12 注入システム



12 メンテナンス用部品 注入システム

インジェクションバルブアセンブリ

品目	部品番号	説明
	0101-1422	インジェクションバルブ
	0100-1852	アイソレーションシール
1	0101-1416	ローターシール PEEK
2	0101-1417	ステータヘッド
3	1535-4857	ステータネジ



カラムオーブン

部品番号	説明
G4280-60040	カラムオーブンアセンブリー式
G4280-60017	ヒータードアアセンブリ



12 メンテナンス用部品 検出器

検出器

可変波長検出器 (VWD)

標準フロー セル 10 mm / 14 µL

品目	部品番号	説明
	G1314-60086	標準フローセル、10 mm、14 μL、40 bar
	5062-8522	キャピラリカラム - 検出器 PEEK、長さ 600 mm、 内径 0.17 mm、外径 1/16 インチ
	G1314-65061	セルリペアキット、内容:2x ガスケット #1、 2x ガスケット #2、2x ウィンドウクォーツ
1	G1314-65062	セルネジキット
2	79853-29100	コニカルスプリングキット、10 個入
3	G1314-65066	#2 リングキット(IN、内径 1 mm の小穴)、PEEK、 数量 = 2
4	G1314-65064	#2 ガスケット(IN、内径 1 mm の小穴)、 カプトン、数量 = 10
5	79853-68742	ウィンドウクォーツキット、2 個入
6	G1314-65063	#1 ガスケットキット (OUT、内径 2.4 mm の大穴)、 カプトン、数量 = 2
7	G1314-65065	#1 リングキット(OUT、内径 2.4 mm の大穴)、 PEEK、数量 = 2



図 74 標準フロー セル 10 mm / 14 µL

12 メンテナンス用部品 検出器

検出器ランプ

部品番号 説明

G1314-60100 重水素ランプ

光学ユニットとファンアセンブリ

部品番号 説明

- G1314-60061 光学ユニットアセンブリー式
- G4280-80004 ファン
- G1314-60114 ヒーターアセンブリ
- G1314-67000 ヒーターインタフェースボードキット(リベット 4 本を含む)
- G1314-65802 VWD 温度センサーボード



注記

光学ユニットの修理には、専門的な知識が必要です。

ダイオードアレイ検出器 (DAD)

標準フロー セル

品目	部品番号	説明
	G1315-60022	標準フローセル、10 mm、13 μL、120 bar(12 MPa)
1	79883-22402	ウィンドウネジ
2	5062-8553	スプリングワッシャ
3	79883-28801	圧縮ワッシャ
4	79883-22301	ウィンドウホルダ
5	1000-0488	クォーツウィンドウ
6	G1315-68711	ガスケット背面 (PTFE)、2.3 mm 穴、 出口側(12 個/pk)
7	G1315-68710	ガスケット前面(PTFE)、1.3 mm 穴、 注入口側(12 個/pk)
8		ウィンドウアセンブリ(ウィンドウネジ、スプリング ワッシャ、圧縮ワッシャ、ウィンドウホルダ、クォー ツウィンドウを含む)
	G1315-87331	キャピラリ IN (0.17 mm、590 mm lg) 熱交換器を含む
10	G1315-87302	キャピラリ OUT (0.17 mm、200 mm)
11	G1315-84910	クランプユニット
	0515-1056	ネジ、M 2.5、4 mm 長 セル本体 / クランプ用
	5022-2184	ユニオン ZDV
	G1315-68712	セル修理キット(標準) ウィンドウネジキットを含む、4 mm 六角レンチおよ びシールキット
	79883-68703	ウィンドウネジキット、クォーツウィンドウ 2 個、 圧縮ワッシャ 2 個、ウィンドウホルダ 2 個、 ウィンドウネジ 2 個、ワッシャ 10 個を含む)



図 75 標準フローセル部品

注記

ガスケット#6と#7は、口径が異なります。



12 メンテナンス用部品 検出器

検出器ランプ

部品番号 説明

- 2140-0820 長寿命重水素ランプ「C」(黒色カバーおよび ID タグ付き)
- G1103-60001 タングステンランプ



この章では、LC システムのアップグレード用の情報を説明します。



13 Agilent 1220 Infinity LC のアップグレード オーブンのアップグレード

オーブンのアップグレード

必要な部品:部品番号説明G4297A1220 Infinity オーブンキット

必要なソフトウェ Lab Advisor ソフトウェア

ア:

- 1 機器をのスイッチを切ります。
- 2 下部フロント カバーを外します。
- 3 カラムを取り外します。
- 4 リークチューブを外します。
- 5 カラムトレイの両側のぎざぎざの部分を内側に押して、カラムトレイを 取り外します。
- **6** オーブンアップグレードキットを開梱し、2 つの部品に分けます。
- 7 オーブンをカラムトレイにカチッとはめ込みます。

オーブンへの電気的接続が自動的に行われます。

- 8 リークチューブを取り付けます。
- 9 オーブンの断熱材を、オーブンの断熱材サポートの切り抜きが下になる ようにして、下部フロント カバーにカチッとはめ込みます。
- 10 カラムを取り付け、キャピラリを再接続します。
- 11 下部前面カバーを取り付けます。
- 12 Lab Advisor ソフトウェアを開始し、機器を接続します。機器コント ロールを開いた後、任意のサブモジュールについて変換を開き、オーブ ン追加をクリックします。ハードウェアのコンフィグレーションの変更 は、次の電源のオン / オフで終了します。
- 13 機器の電源のオン / オフ後、クロマトデータシステムを開始します。自動構成を使用し、これで機器は新しいメインアセンブリ番号下にレジス タされます。



1220 Infinity LC

14 ケーブルの識別

ケーブル概要 394 アナログケーブル 396 リモートケーブル 398 BCD ケーブル 402 CAN/LAN ケーブル 405 Agilent 1200 モジュールから PC 406

この章では、1200 シリーズの HPLC モジュールに使用されるケーブ ルについて説明します。



393

14 ケーブルの識別 ケーブル概要

ケーブル概要

注記

安全規準または EMC 規格のコンプライアンスと正しい動作を確実にするため に、Agilent Technologies 製以外のケーブルは使用しないでください。

アナログケーブル

部品番号	説明
35900-60750	Agilent モジュールから 3394/6 インテグレータ
35900-60750	Agilent 35900A A/D コンバータ
01046-60105	アナログケーブル (BNC から汎用、スペードラグ)

リモートケーブル

部品番号 説明 03394-60600 Agilent モジュールから 3396A シリーズ I インテグレータ 3396 シリーズ II/3395A インテグレータについては、『「リ モートケーブル」 398 ページ』 セクションの詳細を参照して ください。 03396-61010 Agilent モジュールから 3396 シリーズ III/3395B インテ グレータ 5061-3378 リモートケーブル 01046-60201 Agilent モジュールから汎用

BCD ケーブル

部品番号 説明

- 03396-60560 Agilent モジュールから 3396 インテグレータ
- G1351-81600 Agilent モジュールから汎用

CAN ケーブル

部品番号	説明				
5181-1516	CAN ケーブル、	Agilent	モジュール間、	0.	5 m
5181-1519	CAN ケーブル、	Agilent	モジュール間、	1	m

LAN ケーブル

部品番号 説明

5023-0203	クロスオーバーネットワークケーブル、シールド付き、 3 m(ポイントツーポイント接続用)
5023-0202	ツイストペアネットワークケーブル、シールド付き、7 m (ポイントツーポイント接続用)

RS-232 ケーブル

部品番号 説明

G1530-60600 RS-232 ケーブル、2 m

RS232-61601
 RS-232 ケーブル、2.5 m
 機器のPC接続用、9-to-9 ピン(メス)このケーブルのピン
 アウトは特殊で、プリンタやプロッタの接続はできません。
 このケーブルは、書き込みをピン 1-1、2-3、3-2、4-6、
 5-5、6-4、7-8、8-7、9-9 で行う、フルハンドシェークの
 「ヌルモデムケーブル」ともいいます。

5181-1561 RS-232 ケーブル、8 m



アナログケーブル



アナログケーブルの一端は、Agilent モジュールに接続できる BNC コネク タになっています。もう一端は、接続する機器によって異なります。

Agilent モジュールから 3394/6 インテグレータ

部品番号 35900-60750	ピン 3394/6	ピン Agilent モジュール	シグナル名
	1		未接続
	2	シールド	アナログ -
	3	センタ	アナログ +
Agilent モジュールから BNC コネクタ

部品番号 8120-1840	ピン BNC	ピン Agilent モジュール	シグナル名
	シールド	シールド	アナログ -
	センタ	センタ	アナログ +

Agilent モジュールから汎用への接続

部品番号 01046-60105	ピン	ピン Agilent モジュール	シグナル名
	1		未接続
-E	2	黒	アナログ -
	3	赤	アナログ +
- ZE			



リモートケーブル



このタイプのケーブルの一端は、Agilent モジュールに接続できる APG (Analytical Products Group) リモートコネクタになっています。もう一端は、接続する機器によって異なります。

Agilent モジュールから 3396A インテグレータ

部品番号 03394-60600	ピン 3396A	ピン Agilent モジュール	シグナル名	アク ティブ (TTL)
	9	1 - 白	デジタルグ ランド	
	NC	2 - 茶	プレラン	低
	3	3 - 灰	[スタート]	低
	NC	4 - 青	シャットダ ウン	低
	NC	5 - ピンク	未接続	
	NC	6 - 黄	電源オン	青
	5,14	7 - 赤	レディ	青
	1	8 - 緑	ストップ	低
	NC	9 - 黒	スタートリ クエスト	低
	13, 15		未接続	

Agilent モジュールから 3396 シリーズ 11/3395A インテグレー タ

ケーブル Agilent モジュールから 3396A シリーズ I インテグレータ (03394-60600) のインテグレータ側のピン #5 を切断して使用します。切 断しないで使用すると、インテグレータは START; not ready を印字しま す。

Agilent モジュールから 3396 シリーズ 111/3395B インテグレー タ

部品番号 03396-61010	ピン 33XX	ピン Agilent モジュール	シグナル名	アク ティブ (TTL)
	9	1 - 白	デジタルグ ランド	
	NC	2 - 茶	プレラン	低
	3	3 - 灰	[スタート]	低
	NC	4 - 青	シャットダ ウン	低
	NC	5 - ピンク	未接続	
	NC	6-黄	電源オン	青
	14	7 - 赤	レディ	青
	4	8 - 緑	ストップ	低
	NC	9 - 黒	スタートリ クエスト	低
	13, 15		未接続	

部品番号 5061-3378	ピン 35900 A/D	ピン Agilent モジュール	シグナル名	アク ティブ (TTL)
	1 - 白	1 - 白	デジタルグ ランド	
0 50	2 - 茶	2 - 茶	プレラン	低
	3 - 灰	3 - 灰	[スタート]	低
0	4 - 青	4 - 青	シャットダ ウン	低
	5 - ピンク	5 - ピンク	未接続	
	6 - 黄	6 - 黄	電源オン	青
	7 - 赤	7 - 赤	レディ	青
	8 - 緑	8 - 緑	ストップ	低
	9 - 黒	9 - 黒	スタートリ クエスト	低

Agilent モジュールから Agilent 35900 A/D コンバータ

部品番号 01046-60201	ワイアの色	ピン Agilent モジュール	シグナル名	アク ティブ (TTL)
	É	1	デジタルグ ランド	
	茶	2	プレラン	低
	灰	3	[スタート]	低
	青	4	シャットダ ウン	低
	ピンク	5	未接続	
	黄	6	電源オン	高
	赤	7	レディ	高
	緑	8	ストップ	低
	黒	9	スタートリ クエスト	低

Agilent モジュールから汎用への接続



BCD ケーブル



BCD ケーブルの一端は、Agilent モジュールに接続できる 15 ピンの BCD コネクタになっています。もう一端は、接続する装置によって異なります。

部品番号 G1351-81600	ワイアの色	ピン Agilent モジュール	シグナル名	BCD の 桁
	緑	1	BCD 5	20
19	柴	2	BCD 7	80
	青	3	BCD 6	40
	黄	4	BCD 4	10
	黒	5	BCD 0	1
	オレンジ色	6	BCD 3	8
	赤	7	BCD 2	4
	茶	8	BCD 1	2
	灰	9	デジタルグ ランド	灰
	灰 / ピンク	10	BCD 11	800
	赤 / 青	11	BCD 10	400
	白/緑	12	BCD 9	200
	茶/緑	13	BCD 8	100
	未接続	14		
	未接続	15	+ 5 V	低

Agilent モジュールから汎用への接続

部品番号 03396-60560	ピン 3396	ピン Agilent モジュール	シグナル名	BCD の 桁
	1	1	BCD 5	20
80,15	2	2	BCD 7	80
	3	3	BCD 6	40
	4	4	BCD 4	10
	5	5	BCD 0	1
	6	6	BCD 3	8
	7	7	BCD 2	4
	8	8	BCD 1	2
	9	9	デジタルグ ランド	
	NC	15	+ 5 V	低

Agilent モジュールから 3396 インテグレータ

CAN/LAN ケーブル



CAN/LAN ケーブルの両端は、Agilent モジュールの CAN または LAN コネ クタに接続できるモジュラプラグになっています。

CAN ケーブル

部品番号	説明
5181-1516	CAN ケーブル、Agilent モジュール間、0.5 m
5181-1519	CAN ケーブル、Agilent モジュール間、1 m

LAN ケーブル

部品番号	説明
5023-0203	クロスオーバーネットワークケーブル、シールド付き、 3 m(ポイントツーポイント接続用)
5023-0202	ツイストペアネットワークケーブル、シールド付き、7 m (ポイントツーポイント接続用)

14 ケーブルの識別

Agilent 1200 モジュールから PC

Agilent 1200 モジュールから PC

部品番号 説明

G1530-60600 RS-232 ケーブル、2 m

- RS232-61601
 RS-232 ケーブル、2.5 m
 機器の PC 接続用、9-to-9 ピン (メス) このケーブルのピン
 アウトは特殊で、プリンタやプロッタの接続はできません。
 このケーブルは、書き込みをピン 1-1、2-3、3-2、4-6、
 5-5、6-4、7-8、8-7、9-9 で行う、フルハンドシェークの
 「ヌルモデムケーブル」ともいいます。
- 5181-1561 RS-232 ケーブル、8 m



1220 Infinity LC

15 付録

安全に関する一般的な情報 408 溶媒情報 411 無線干渉 413 紫外線放射 414 騒音レベル 415 廃電気電子機器 (WEEE) 指令 (2002/96/EC) 416 ホルミウムフィルタの波長の証明書 417 Agilent Technologies のインターネットサービス 418



安全に関する一般的な情報

安全に関する一般的な情報

安全に関する一般的な情報

以下の安全に関する一般的な注意事項は、本機器の操作、サービス、および修理のすべての段階で遵守するようにしてください。以下の注意事項またはこのマニュアルの他の箇所に記載されている警告に従わないと、本機器の設計、製造、および意図された使用法に関する安全基準に違反することになります。使用者側による遵守事項からのかかる逸脱に起因する問題について Agilent は免責とさせて頂きます。

警告 装置の正しい使用法を確保してください。 機器により提供される保護が正常に機能しない可能性があります。

→この機器のオペレーターは、本マニュアルで指定した方法で機器 を使用することをお勧めします。

安全規格

本製品は、国際安全基準に従って製造および試験された、安全クラス I 装置(アース端子付き)です。

操作

電源を投入する前に、設置方法が本書の説明に合っているかどうか確認してください。さらに、次の注意を守ってください。

操作中に装置のカバーを取り外さないでください。装置のスイッチを ON に する前に、すべての保護接地端子、延長コード、自動変圧器、および本装 置に接続されている周辺機器を、接地コネクタを介して保護接地に接続し てください。保護接地がどこかで途切れていると、感電によって人体に重 大な危害を及ぼすことがあります。保護が正常に機能していないと思われ る場合は、装置のスイッチを OFF にして、装置の操作を中止してください。

安全に関する一般的な情報

ヒューズを交換する際は、必ず指定したタイプ(普通溶断、タイムラグなど)と定格電流のヒューズだけを使用してください。修理したヒューズを 使用したり、ヒューズホルダを短絡させたりしてはなりません。

本書で説明した調整作業には、装置に電源を入れた状態で、保護カバーを 取り外して行うものがあります。その際に、危険な箇所に触れると、感電 事故を起こす可能性があります。

機器に電圧をかけた状態で、カバーを開いて調整、メンテナンス、および 修理を行うことは、できるだけ避けてください。どうしても必要な場合は、 経験のある担当者が感電に十分に注意して実行するようにしてください。 内部サービスまたは調整を行う際は、必ず応急手当てと蘇生術ができる人 を同席させてください。メンテナンスを行うときは、必ず装置の電源を 切って、電源プラグを抜いてください。

本装置は、可燃性ガスや有毒ガスが存在する環境で操作してはなりません。このような環境で電気装置を操作すると、引火や爆発の危険があります。

本装置に代替部品を取り付けたり、本装置を許可なく改造してはなりません。

本装置を電源から切り離しても、装置内のコンデンサはまだ充電されてい る可能性があります。本装置内には、人体に重大な危害を及ぼす高電圧が 存在します。本装置の取り扱い、テスト、および調整の際は十分に注意し てください。

特に、有毒または有害な溶媒を使用する場合は、試薬メーカーによる物質 の取り扱いおよび安全データシートに記載された安全手順(保護眼鏡、安 全手袋、および防護衣の着用など)に従ってください。

安全に関する一般的な情報

安全記号

表 48 安全記号

記号	説明
\wedge	危害のリスクを保護するために、そして装置を損傷から 守るために、ユーザーが取扱説明書を参照する必要があ る場合、装置にこの記号が付けられます。
4	危険電圧を示します。
	アース (保護接地) 端子を示します。
	本製品に使用されている重水素ランプの光を直接目で見 ると、目をいためる危険があることを示しています。
<u>s</u>	表面が高温の場合に、この記号が装置に付けられます。 加熱されている場合はユーザーはその場所を触れないで ください。

警告は、

警告

注意

人身事故または死に至る状況を警告します。

→ 指示された条件を十分に理解してそれらの条件を満たしてから、 その先に進んでください。

注意

データ損失や機器の損傷を引き起こす状況を警告します。

→ 指示された条件を十分に理解してそれらの条件を満たしてから、その先に進んでください。

溶媒情報

溶媒を使用するときは、次の注意に従ってください。

フローセル

アルカリ性溶液 (pH > 9.5) は、クォーツを侵し、フローセルの光学特性を劣化させるため、使用を避けてください。

バッファ溶液が結晶化しないようにします。これにより、フローセルの詰 まりや損傷を引き起こします。

フローセルを 5°C より低い温度で輸送する場合は、必ずセルにアルコールを満たします。

フローセル内の水性溶媒は、藻を増やすことがあります。そのため、フ ローセル内に水性溶媒を残さないでください。数%の有機溶媒(たとえ ば、アセトニトリルまたはメタノール~5%)を添加します。

溶媒

褐色の溶媒ボトルを使用すると藻の発生を避けることができます。

微粒子による配管の詰まりを避けるために、溶媒は必ずろ過します。また、 次の鉄腐食性溶媒の使用は避けてください。

- ハロゲン化アルカリ化合物およびその酸溶液(ヨウ化リチウム、塩化カ リウムなど)。
- 硝酸、硫酸などの高濃度の無機酸(特に高温の場合)。(可能な限り腐食 性の弱いリン酸またはリン酸緩衝液に変更して下さい)。
- 以下に示すラジカルまたは酸、あるいはその両方を発生するハロゲン化 溶媒または混合液。

 $2CHC1_3 + 0_2 \rightarrow 2COC1_2 + 2HC1$

乾燥プロセスによって安定剤のアルコールが除去された場合、通常はス テンレスを触媒として、乾燥クロロホルムでこの反応が急速に発生しま す。

- THF、ジオキサン、ジイソプロピルエーテルなどのクロマトグラフィグレードのエーテルは過酸化物を含む可能性があります。このようなエーテルは、過酸化物を吸収する乾性アルミニウム酸化物でろ過してください。
- 有機溶媒中の有機酸溶液(酢酸、ギ酸など)。たとえば、酢酸1%のメタノール溶液は鋼鉄を侵します。
- ・ 強力なキレート試薬(EDTA、エチレンジアミン 4 酢酸など)を含む溶液。
- ・四塩化炭素と 2- プロパノールまたは THF の混合溶液。

無線干渉

無線干渉に対して最適な保護を行うために、アジレントが提供するケーブルは選別されています。すべてのケーブルが安全性または EMC 規格に準拠しています。

テストと測定

選別していないケーブルを用いてテスト機器と測定機器を操作したり、確定していない設定での測定に使用する場合、無線干渉が制限する運転条件がまだ許容範囲内であることをユーザーが確認する必要があります。

紫外線放射

紫外線放射

本製品による紫外線放射 (200 - 315 nm) の制限値は、オペレータや点検 作業者の防護されていない皮膚または目に対する暴露について、American Conference of Governmental Industrial Hygienists (米国産業衛生専門 家会議) により既定される以下の TLV (限界値) を遵守するものです。

表 49 UV 放射リミット

暴露時間 / 日	有効放射照度
8 h	0.1 μ W/cm ²
10 min	5.0 μ W/cm ²

通常の放射量は、これらのリミットを大幅に下回ります。

表 50 UV 放射標準値

ポジション	有効放射照度
据付済みランプ、距離 50 cm	平均 0.016 µW/cm ²
据付済みランプ、距離 50 cm	最大 0.14 µW/cm ²

騒音レベル

製造業者による宣言

本製品は、ドイツ騒音条例(1991 年 1 月 18 日)の条件に適合しています。

本製品の音圧レベル(オペレータの位置)は、70 dB 未満です。

- 音圧 L_p < 70 dB (A)
- オペレータの位置
- 通常動作時
- ISO 7779:1988/EN 27779/1991 (タイプテスト) に準拠

廃電気電子機器(WEEE)指令(2002/96/EC)

廃電気電子機器(WEEE)指令(2002/96/EC)

要約

2003 年 2 月 13 日に欧州委員会が可決した、廃電気電子機器(WEEE) 指令(2002/96/EC) は、すべての電気および電子機器に関する生産者責任を2005 年 8 月 13 日から導入するというものです。

注記 本製品は、WEEE 指令(2002/96/EC)に準拠しており、要件を記しています。 貼り付けられたラベルには、この電気 / 電子機器を家庭用廃棄物として廃棄し てはならないことが表示されています。

製品カテゴリ:

WEEE 指令付録 I の機器の種類を参照して、本製品は「モニタリングおよび 制御装置」製品と分類されます。



注記

家庭用廃棄物として捨ててはいけません

不必要な製品を返品するには、最寄りのアジレント営業所にお問い合わせいただくか、詳細について www.agilent.com をご覧ください。

ホルミウムフィルタの波長の証明書

ホルミウムフィルタの波長の証明書

Declaration of Conformity					
We herewith inform you that the					
Holmium Oxide Glass Filter					
used in Agilents absorbance detectors listed in the table below meets the requirements of National Institute of Standards and Technology (NIST) to be applied as certified wavelength standard.					
According to the publication of NIST in J. Res. Natl. Inst. Stand. Technol. 112, 303-306 (2007) the holmium oxide glass filters are inherently stable with respect to the wavelength scale and need no recertification. The expanded uncertainty of the certified wavelength values is 0.2 nm.					
Agilent Technologies guarantees, as required by NIST, that the material of the filters is holmium oxide glass representing the inherently existent holmium oxide absorption bands.					
Test wavelengths:					
Product Number	Series	Measured	Wavelength	Optical	
70000 4	1000	Wavelength	* Accuracy	Bandwidth	
79883A	1090	361.0 nm	+/- 1 nm	2 nm	
/9854A	1050	418.9 nm			
G1300A	1000	435.7 mm			
G1315A, G1365A	1100 / 1200 / 1260	550.7 1111			
G1500A G7100A	CE	+			
70853C	1050	360.8nm	+/- 2 nm	6.000	
190550	1050	418 5nm	·/- 2 mm	·	
		536.4nm			
G1314A/B/C	1100 / 1200 / 1260	360.8nm	+/- 1 nm	6 nm	
G1314D/E/F		418.5nm			
G4286,, 90A/B/C	1120 / 1220	536.4nm			
*) The variation in Measured Wavelength depends on the different Optical Bandwidth.					
May 19, 2010					
/	(Date)				
Thomas for					
(R&D Manager) (Quality Manager)					
P/N 89550-90501	Revision: H Effective by: May 19, 2010		Agilent Technologies		

Agilent Technologies のインターネットサービス

Agilent Technologies のインターネットサービス

製品およびサービスの最新情報を知るには、アジレントのウェブサイトに アクセスしてください。

http://www.agilent.com

Products/Chemical Analysis を選択してください。

索引

A

ASTM 環境条件 18 基準と条件 26 基準 29 AUTO モード 94

В

BCD ケーブル 402 Beer-Lambert (法則) 141 Bootp サービス 停止 76 インストール 69 再起動 77 設定 76 Bootp を使用した自動コン フィグレーション 67 Bootp 58 および保存 59 自動コンフィグレーショ ン 67 初期化モード 58 設定の永久保存 78 デフォルトの使用 60 保存されたパラメータの使 用 60

С

CAN 通信消失 217 CAN ケーブル 405

D

D/A コンバータ 202 DAC Agilent Lab 202 Advisor インスタント パイロッ 203 ト DHCP セットアップ 64 一般情報 62 DSP が実行されていませ ん 274

E

Early maintenance feedback 20 EE 2060 227 EMF カウンタ、ポンプ 12 ポンプヘッド 12

G

GLP 機能 20

L

Lab Advisor 279 LAN Bootp および保存 59 Bootp を使用した自動コン フィグレーション 67 Bootp 58

TCP/IP パラメータコン フィグレーション 56 telnet を使用したマニュ アルコンフィグレーショ ン 80 ケーブル 405 初期化モード選択 58 設定の永久保存 78 デフォルトの使用 60 保存されたパラメータの使 用 60 マニュアルコンフィグレー ション 79 リンクコンフィグレーショ ン選択 66 LMD 20

Μ

MAC アドレス 確認 73 MCGV ヒューズ 234 message heater failed 252

Ρ

pH 範囲 21, 22 PM 279 PTFE フリット 290

R

RFID (無線認識) タ グ 122 RS-232C

1220 Infinity LC

索引

S SSV 10, 11 T TCP/IP パラメータコンフィグ レーション 56 telnet コンフィグレーショ ン 80

ケーブル 406

U

URL 418 UV ランプオン 179 UV 放射 414

۷

VWD EMF カウンタ 14

W

WEEE 指令 416

Х

X-軸 109

Z

Z- 軸 109

あ

アーム移動 161 アームを上げる 164 アームを下げる 164 アームをホームへ移 動 161

アクロマート ソース レンズ 123 アレンジト・テクノロ ジー 418 アジレントメンテナン ス 279 アダプタ 286 アナリティカルヘッ ド 107 アナログ ケーブル 396 アライメント グリッパ 167 アライメントティーチン グ 164 アルファおよびベータ ライ ン 199 アレイ 124 圧力 86 圧縮率補正 21, 22, 92 圧力、動作範囲 21. 22 圧力が下限値より低 い 226 圧力が上限値を超過 225 圧カシグナルが読み取れな い 227 圧力測定値が読み取れな い 223 圧力の脈動 21. 22. 92. 94 圧力プロファイル 149. 150 安全クラス 丨 408 安全機能 システム 20 安全 規格 19

記号 410 一般的な情報 408 暗電流 171 テスト 185

い

インジェクションシー ル 304 インジェクションシール vespel 99 インジェクションルール tefzel 99 インターネット 418 インデックスが見つからな い 222 インデックス調整 220 インデックスリミッ F 221 移送機構 102 移動相と接触する部品の材 質 90. 91 情報 キュベットホルダ 335 一般エラーメッセー ジ 209

え

エラー メッセージ DSP が実行されていません 274 uv 点灯失敗 269 uv ヒータ電流 270 uv ランプ電圧 256 uv ランプ電流 255 グレーティングがありません 262

ダイオード電流リー 268 ク 波長キャリブレーション失 敗 271 波長リキャリブレーション 消失 273 フィルタがありませ 260 ん ポンプ コンフィグレー ション 227 エラーメッセージ ADC ハードウェアエ ラー 257 lost CAN partner 217 MCGV ヒューズ 234 アーム移動失敗 237 圧力が上限値より低 い 226 圧力が上限値を超 過 225 圧カシグナルが読み取れな い 227 圧力測定値が読み取れな い 223 安全フラップ不良 248 インデックスが見つからな い 222 インデックス調整 220 インデックスリミッ F 221 エンコーダが見つからな い 219 温度上限を超過 231 キャリブレーション失 敗 258 空気吸入口のセンサからの 不正な温度値 254 グリッパ内にバイアルあ り 251

グレーティング / フィルタ モーター故障 261 サーボ再起動の失 229 敗 シャットダウン 216 初期化失敗 223. 239 ストローク長 230 洗浄用バイアルが見つから ない 244 タイムアウト 218 タイムアウト待ち 235 ニードルアップの失 敗 247 ニードルダウンの失 246 敗 バイアル初期化 240 バイアルなし 243 バイパスへのバルブ切り替 え失敗 249 波長チェック失敗 266 バルブ故障 233 範囲外の温度 232 ヒーター故障 252 ヒーター出力限界 253 ヒーター電流の不 足 263 ファンアセンブリのセンサ からの不正値 253 ファン動作不良 211 フィルタチェック失 敗 259 補正センサーオープ 209 ン 補正センサーショー 210 ト ホルミウムオキサイドテス ト失敗 264 ポンプヘッドが見つからな 228 い

メインパスへのバルブ切り 替え失敗 250 モーターの障害 245 モータドライブ電 カ 224 溶媒ゼロカウンタ 236 ランプ点火不良 265 リークセンサーオープ ン 213 リークセンサーショー ト 214 212 リーク リモートタイムアウ 215 ト 計量ピストンのホームポジ ションへの移動失 敗 242 無効なバイアルポジショ 241 ン エンコーダが見つからな い 219 液量メータ 12 お オート サンプラ EMF カウンタ 13 簡単な修理 307 概要 102 オートサンプラ オートサンプラアセンブリ 部品 308 オートサンプラのクリーニン グ 309 オート サンプラの取り付 け サンプル トレイ 110 オーブンキャリブレーショ ン 169

索引

オーブンテスト 168 温度 23 温度上限を超過 231 温度センサー 212

か

カウンタ インジェクションバル ブ 13 総液量 12 ニードル動作 13 シール磨耗 12 オート サンプラ 13 検出器 14 カバーなし再スター F 228 概要 光学ユニットの部 品 123 可変磁気抵抗モータ 88 可変ストローク容量 94 124 可変入射スリット カラムオーブン 118 環境 18 緩衝液 300 簡単な修理 オート サンプラ 307

き

キャップ 111 キャリブレーション オーブン 169 キュベットホルダ 335 基準条件 ASTM 29 吸引 162 吸光度 Beer-Lambert 141 強度テスト 173, 187 凝縮 18

<

グラジエント形成 21. 22 グラジエントバルブ (DCGV) 300 グリッパ 109 交換 161 グリッパアーム 修理 325 グリッパフィンガー 109 グリッパベリフィケーショ ン 167 グリッパを閉じる 165 グリッパを開く 165 グレーティング 124. 124 グレーティングモー ター 176

け

ケーブル BCD 402. 395 CAN 405. 395 LAN 405. 395 RS-232 406. 395 アナログ 396. 394 概要 394 リモート 398. 394 検出タイプ 25 検出器 EMF カウンタ 14 121 特徴

検出 化合物クラス 135 E コントロール システム 20 コンフィグレーション 10 高圧異常チェック 157 交換 インジェクションシー 304 ル グリッパ 161 デュアル チャンネル グラ ジェントバルブ (DCGV) 300 ニードル 160 パージ バルブ 290 パージバルブフリッ 290 ۲-パッシブインレットバル ブ 286 ピストン 160 梱包明細リスト 32

さ

サーボ再起動の失敗 229 サファイア ピストン 88. 88 サンプルおよびリファレンス 波長 128 サンプルスキャン 179 サンプル トレイ バイアルポジションの番号 付け 110 サンプルトレイ 110 サンプルの注入 98 サンプル量 99

サンプル ループ 98 サンプリング動作 103 サンプリングユニッ F 106 最適化 サンプルおよびリファレン ス波長 128 スペクトル取り込 み 133 スリット幅 131 選択性の 134 ピーク幅 126 酸化ホルミウム フィルタ 123

L

シーター軸 109 シグナルプロット 148 システム圧力上昇 360 システムコンフィグレーショ ン 10 シャットダウン 216 シール摩耗カウンタ 12 湿度 19 周波数範囲 19 修理方法 インジェクションシー ル 304 仕様 23. 29 性能 25 ダイオード幅 27 直線範囲 27 ノイズおよび直線 29 性 ノイズおよびドリフト (ASTM) 27 波長真度 27

波長範囲 27 物理的 19. 19 フローセル 28 プログラム式スリット 幅 27 初期化失敗 223 初期化モード選択 58 初期化 ポンプ 90 診断テーブル 148 修理 簡単な修理の概要 327 機器のクリーニン グ 339 ニードルアセンブ IJ 310 ファームウェアの交 換 361 メタリングシール 321 メタリングプラン ジャ 321 リーク処理システムの交 358 換 リークの補正 357 ロータ シール 317 手動コントロール 161 使用 キュベットホルダ 335 使用温度 19 使用高度 19 使用周囲温度 19 消費電力 19 重量 19 情報 148 UV 放射について 414 溶媒について 411 修理

ニードルシートアセンブ リ 315

す

スキャン 1	79	
ステッパモータ	107	
ステータス情報	t 148	
ステータ 1	08	
ステップコマン	ド 161	
ストローク長	230	
ストローク容量	88, 94	
スペクトルスキ	・ヤン 179	
スペクトル		
取り込み	133	
スリット幅	27, 131	
推奨 pH 範囲	21, 22	
水平部、リーク	テス	
F 154		
据え付けチェッ	ク 147	
据付		
1/H I J		
梱包明細リン	スト 32	

せ

セル 170 静電気放電(ESD) 281 精度 99 性能仕様 23 性能 仕様 25 設置スペース 17 設置について 16 環境 18 設置スペース 17 電源コード 17 電源について 16

索引

設置

設置について 設定可能な流量範囲 21, 22 設定の永久保存 31 説明 98 選択性の最適化 134

そ

測光精度 142 組成精度 21, 22 組成範囲 21, 22

た

タイムアウト 218 タイムアウト待ち 235 ダイオード アレイ 125, 124 幅 27

ち

注入ロバルブ 286 注入バルブ 102, 108 注入 98, 100 注入動作 104 直線性 25, 26 仕様 29 直線範囲 27

つ

通信 20 詰まり 157, 224

て

テスト機能 D/A コンバータ 202

DAC 202 テストクロマトグラ ム 178, 197 テスト 暗電流 185 強度 187 テストクロマトグラ ム 197 波長キャリブレーショ ン 177 フィルタ 182 ホルミウムオキサイ ド 190 ディレイ ボリューム 21. 90. 91 データのエクスポー F 179 データ評価 20 デガッサ 95 デュアル チャンネル グラジ エントバルブ 300 デュアルプランジャ直列型設 計 86 定量バルブ 高速 86 電圧範囲 19 電源周波数 19 電源 コード 17 留意事項 16

لح

トラブルシューティング エラーメッセージ 208 トランスポート アセンブ リ 109 ドリフト 25 ドリフト (ASTM) およびノイ ズ 27 動作可能圧力範囲 21, 22

に

ニードル 101 交換 160 ニードルアップ 162. 162 ニードルタイプ 101 ニードルドライブ 106. 107 ニードルをサンプルに挿 入 162 ニードルをシートに挿 入 162 入射スリット 124 入力電圧 19

の

ノイズおよび直線性 仕様 29 ノイズおよびドリフト (ASTM) 27

は

ハーフトレイ 110
ハイドロリックシステム 21, 22
ハイドロリックパス 86
バイアル内容物温度 23
バイアルの番号付け 110, 110
バイアル 102, 111
バイアルラック 102

1220 Infinity LC

バイアルをシートに移 162 動 バイアルをトレイ戻 す 162 バイパス 104 バッファアプリケーショ ン 95 バルブ故障 233 バルブのメインパス 163 バルブバイバス 162 バルブ 定量 86 バルブフリット 290 バンド幅 6.5 nm 25 パージ バルブ 290 パージバルブフリット 95 パージポンプ 159 廃棄物 電気電子機器 416 廃電子機器 416 波長ベリフィケーションおよ びリキャリブレーショ ン 199 波長リキャリブレーション消 失 273 波長 キャリブレーショ 177 ン 直度 25. 27 範囲 190 ~ 600 25 nm 範囲 27 ベリフィケーションおよび リキャリブレーショ ン 199 範囲外の温度 232

ひ ピーク幅(レスポンス タイ ム) 126 ピストンチャンバ 86 ピストン 88, 95 交換 160

ふ

ファームウェア 148 アップグレード / ダウング レード 361. 361 更新 361 更新 361 ファン動作不良 211 フィルタ / グレーティングテ スト 176 フィルタ テスト 182 フィルタモーター 176 フローセル 140 タイプとデータ 25 フローセルの補正ファク 142 タ フローセル 123. 170 サポートウィンド ゥ 123 仕様 28 384 標準(部品) 補正ファクタ 142 ブランクスキャン 179 プランジャホーム 162 プログラム式スリット 幅 27 物理的仕様 19. 19 部品の識別 ケーブル 393 分光器 123, 124

nm 当りのダイオー ド 124

^

ベータおよびアルファ ライ ン 199

ほ

ホルミウムオキサイド テスト 190 ボールスクリュードライ ブ 88 $\pi - 1 \sim 106$ ポンプ コンフィグレーショ ン 227 ポンプの異常 359 ポンプピストン 95 ポンプヘッドが見つからな い 228 ポンプヘッドの再設 299 置 ポンプヘッド 再設置 299 ポンプ 86 うまく使用するためのヒン 95 ト 86 概要 機能原理 88 ポンプリークテスト 149 保管温度 19 保管高度 19 保管周囲温度 19 補正センサーオープ ン 209 補正センサーショー ト 210

索引

マ

マニュアルコンフィグレー ション LAN 79 マルチ注入オプショ ン 102

む

無線妨害 413

め

メインパス 104 メイク - ビフォア - ブレイ ク 98 メタリングデバイス 106 メッセージ ADC ハードウェアエ ラー 257 uv ヒータ電流 270 uv 点灯失敗 269 uv ランプ電圧 256 uv ランプ電流 255 キャリブレーション失 敗 258 キャリブレーション消 266 失 空気吸入口のセンサからの 不正な温度値 254 グレーティング / フィルタ 261 モーター故障 グレーティングがありませ 262 ん ダイオード電流リー ク 268 波長キャリブレーション失 敗 271 266 波長チェック失敗 ヒーター出力限界 253

ヒーター電流の不 足 263 ファンアセンブリのセンサ からの不正値 253 フィルタがありませ 6 260 フィルタチェック失 259 敗 ホルミウムオキサイドテス ト失敗 264 ランプ点火不良 265 リモートタイムアウ ト 215 メンテナンス機能 309 メンテナンス機能 ステップコマンド 161 メンテナンス キュベットホルダの使 335 用 概要 338 ファームウェアの交 換 361 ファームウェアの交 換 361 328 ランプの交換 メンテンス用部品 標準フロー セル 384

ŧ

モータドライブ電力 224 モジュールオプショ ン 148 モジュール情報 148 モジュールのクリーニン グ 339 藻の情報 411 藻の繁殖 96

よ 溶媒インレットフィル 95 タ 溶媒切り替えバルブ 10. 11 溶媒情報 411 溶媒ゼロカウンタ 236 溶媒送液システム 86 溶媒フィルタ クリーニング 285 チェック 284 詰まり防止 96

6

ランプ強度 173 ランプ 123 タイプ 25

り

リーク テストの実行 152 リークセンサーオープ ン 213 リークセンサーショー ト 214 リークテスト 150 アイソクラティックポン プ 149 リーク電流 171 リーク 212 補正 357 リセット 163 リモート ケーブル 398 流量精度 21. 21. 22. 22 流量範囲 21. 22

流量

不安定 359 流路の詰まり 157 リンクコンフィグレーション 選択 66

れ

レスポンス タイム(ピーク 幅) 126

3

ロータ シール 交換 317 ロード 98, 99 www.agilent.com

本書の内容

本書には Agilent 1220 Compact LC システム を使用、メンテナンス、修理、アップグレー ドする方法に関する情報が記載されていま す。

本書は以下の章で構成されています。

- 概要
- 据付
- ・ Agilent 1220 Infinity LC の説明
- テスト機能とキャリブレーション
- エラー情報
- メインテナンスと修理
- ・ メンテナンスおよび修理用部品
- ・ Agilent 1220 Infinity LC のアップグ レード
- 付録

© Agilent Technologies 2010-2012

Printed in Germany 05/2012



G4280-96016

