



マニュアル





Agilent Technologies



© 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

G4220-96006 Rev. B

エディション

07/2012

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 1290 Infinity バイナリポンプ (G4220A) について説 明します。

1 はじめに

この章では、ポンプの概要、機器の概要、内部コネクタについて説明しま す。

2 設置について

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の情報を示します。

3 モジュールの設置

この章では、所有するシステムで推奨されるスタックのセットアップと、 Agilent 1290 Infinity バイナリポンプの設置についての情報を示します。

4 ポンプの使用

この章では、Agilent 1290 Infinity バイナリポンプの操作パラメータに ついて説明します。

5 性能の最適化

この章では、追加デバイスの性能または使用の最適化のヒントについて説 明します。

6 トラブルシューティングおよび診断

トラブルシューティングと診断機能についての概要

7 エラー情報

この章では、エラーメッセージの意味を解説し、考えられる原因に関する 情報とエラー状態から回復するための推奨方法を示します。

本書の内容

- 7スト機能とキャリブレーション
 この章では、モジュールのテストについて説明します。
- 9 メンテナンス

この章では、Agilent 1290 Infinity バイナリポンプのメンテナンスについて説明します。

10 メンテナンス用部品と器材

この章では、メンテナンス用部品について説明します。

11 ケーブルの識別

本章では、すべてのケーブルに関する情報を記載します。

12 ハードウェア情報

この章では、ハードウェアと電子機器に関してポンプの詳細を説明します。

13 LAN コンフィグレーション

この章では、検出器の Agilent ChemStation PC への接続について説明します。

14 付録

この章では、安全性、法律、ウェブに関する追加情報を記載しています。

1 はじめに 9

機能 10 バイナリポンプの概要 11 ポンプの原理 12

2 設置について 15

設置について 16 物理的仕様 19 性能仕様 20

3 モジュールの設置 23

モジュールの開梱 24
スタック構成の最適化 26
輸送用緩衝材の取り外し 31
ポンプの設置 32
ポンプへの配管 36
シールウォッシュオプションの取り付け 40

4 ポンプの使用 41

バイナリポンプの準備 42 HPLC システムでの藻の繁殖 43 機器コントロール インターフェイスでのポンプの設定 45 ポンプのプライミング 58 溶媒情報 60

5 性能の最適化 67

ディレイボリュームとエクストラカラムボリューム 68 最適ディレイボリュームの設定方法 69 高分解能を達成する方法 71 溶媒キャリブレーションテーブルの使用 74

6 トラブルシューティングおよび診断 77
 モジュールのインジケータとテスト機能の概要 78
 ステータスインジケータ 79
 使用可能なテストとユーザーインタフェース 81
 Agilent Lab Advisor ソフトウェア 82

7 エラー情報 83 エラーメッセージ内容 85 一般エラーメッセージ 86 ポンプのエラーメッセージ 96

8 テスト機能とキャリブレーション 119
 はじめに 120
 システム圧力テスト 121
 ポンプヘッドのリークテスト 124

9 メンテナンス 127

メンテナンスの概要 128 警告と注意 130 メンテナンスの概要 132 モジュールのクリーニング 133 フィッティングとキャピラリの設置 134 シャットオフバルブパネルの交換 135 圧力センサーの交換 137 インレットバルブの交換 139 アウトレットバルブの交換 141 溶媒切り替えバルブ(SSV)の交換 144 構成の変更または Jet Weaver の交換 146 シールウォッシュポンプの交換 149 スタックインレットバルブの取り外し 150 ポンプヘッドの交換 153 ポンプヘッドの分解 159 プライマリポンプヘッドの分解 161 セカンダリポンプヘッドの分解 164 ポンプヘッドの組み立て 167 パージバルブヘッドの交換 177 高圧フィルタアセンブリの部品の交換 180

バルブレールキットの設置 182	
主電源ヒューズの交換 183	
モジュールファームウェアの交換	185
ポンプモジュールの移動の準備	186

10 メンテナンス用部品と器材 191

メンテナンス部品の概要 192 キャピラリ 193 溶媒キャビネットキット 194 シールウォッシュオプション 195 ポンプヘッドアセンブリの部品 196 プライマリポンプヘッドの部品 198 セカンダリポンプヘッドの部品 202 パージバルブ 206 カバー部品 207 リーク部品 208 ヒューズ 208 アクセサリキット 209 その他 210

11 ケーブルの識別 213

ケーブル概要 214 アナログケーブル 216 リモートケーブル 218 BCD ケーブル 222 CAN/LAN ケーブル 225 RS-232 ケーブルキット 226 Agilent 1200 モジュールからプリンタ 227

12 ハードウェア情報 229

ファームウェアについて 230 電気的接続 233 インタフェース 235 8 ビットコンフィグレーション スイッチの設定 243 機器レイアウト 249 EMF (Early Maintenance Feedback) 250 13 LAN コンフィグレーション 251

最初の必要事項 252 TCP/IP パラメータコンフィグレーション 254 設定スイッチ 255 初期化モード選択 257 動的ホストコンフィグレーションプロトコル (DHCP) 262 リンクコンフィグレーション選択 266 Bootp を使用した自動コンフィグレーション 267 マニュアルコンフィグレーション 278 PC およびユーザーインターフェイス ソフトウェアの設定 284

14 付録 287

安全に関する一般的な情報 288 廃電気電子機器(WEEE)指令(2002-96-EC) 291 無線妨害 292 騒音レベル 293 アジレントのウェブサイト 294



この章では、ポンプの概要、機器の概要、内部コネクタについて説 明します。



機能

このバイナリポンプは、高性能を発揮し、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:

- シールウォッシュ機能により、バッファーを使用したアプリケーション においてポンプシールの寿命を延ばすためにオプションのシールウォッシュ機能の搭載が可能。
- ・ メソッドの柔軟性をもたらす溶媒切り替えバルブ。
- Jet Weaver (ミキサー) により最小限のディレイボリュームで優れたミキシング性能を実現。
- 自動パージ機能によりシステムパージを自動的に実施可能。
- 弾性や溶媒容量変更による影響(圧縮率、熱膨張、混合収縮)を補正するための送液サイクルの自動チューニング機能を搭載。
- 最適な流量真度と組成真度が得られる溶媒の選択が可能。
- ・ 送液パフォーマンスを向上する高速排出機能。
- ・ 信頼性の高い 2 ピストン直列設計。
- スムースで確実な動作を実現する 高分解能のピストンコントロール。
 仕様については、『20ページ 表 2』を参照してください。
- 注記 このバイナリポンプは、Agilent 1290 Infinity 液体クロマトグラフとともに 導入されました。

バイナリポンプの概要

Agilent 1290 Infinity バイナリポンプは、1 つのハウジングに内蔵され た 2 台のポンプから構成されます。高圧混合での 2 液グラジエント送液 が可能です。最高の検出器感度を得るために、特に低流量で最高の流量安 定性が必要なアプリケーションに対して、脱気ユニットを装備しています。 ポンプの流路は、グラジエントの待ち時間が最小になるように最適化され ています。このポンプは高分離能 2.1 mm ID カラムによるハイスループッ ト高速グラジエント分析に適しています。0.05 -5 mL/min で耐圧 1200 bar での送液が可能です。溶媒切り替えバルブ により、チャンネル あたりに 2 つの溶媒の中の 1 つの溶媒からバイナリ混合(アイソクラ ティックまたはグラジエント)を行うことが可能です。高濃度緩衝液に使 用するために、アクティブシールウォッシュ(オプション)が利用可能で す。

11

ポンプの原理

バイナリポンプは2 チャンネルのデュアルプランジャ直列型で、溶媒送液 システムに要求されるすべての必須機能を装備しています。2 つのポンプ アセンブリにより、溶媒の計量と高圧側への送液を行います(最大圧力 1200 bar)。

各チャンネルは、各ピストン用の独立したポンプドライブ、ポンプヘッド、 インレットバルブ、アウトレットバルブ、溶媒熱交換器、アウトレット フィルタで構成されています。2 つのチャンネルは、自動パージバルブと Jet Weaver ミキサの低容量混合部に接続され、35 または100 μ1 のボ リュームを、最適な混合性能を得るために下流で追加することができます。

ポンプ圧を監視するためのシステム圧力センサがパージバルブに接続されています。通常は、ディレイボリュームを最小限に抑えるために、混合部の前にあるポンプの B チャンネルに接続されます。



図 1 配管

ユーザーインタフェースでは、最適なフローと混合精度が得られるように、 各チャンネルで使用する溶媒を選択できます。リップルが低い場合は適切 な溶媒を選択する必要はありませんが、特に圧力範囲全体で高い混合精度 を得るためには、補正アルゴリズムで適切な溶媒特性を選択することが非 常に重要です。溶媒の密度は圧力の影響により大きくなり、排出された特 定の溶媒は、圧力の解放時にカラム全体などに広がります。サンプルが検

出器を通過するときにフローのボリュームを適切な値にするには、密度に 関連するフローの不確実性をポンプモジュールで補正する必要があります。

できる限り安定した圧力を常に提供するために、ポンプで継続的に送液サ イクルを調整し、弾性およびデッドボリューム効果を補正しています。こ の機能によってポンプは安定した正確なフローを提供することができ、個 別のキャリブレーション設定が不要になります。コントロールおよび補正 アルゴリズムのもう 1 つの機能としてリーク補正があります。この機能を 使用すると、ポンプ性能に影響を与えずに、プライマリポンプチャンバ(インレットバルブおよびシール)のわずかなリークを補正することができ ます。

ポンプの堅牢性を向上させる目的で、プライマリピストンの送液時間を短 縮する高速排出機能を使用しているため、リークの実質的な影響を大幅に 軽減することができます。さらに、各ポンプチャンネルのバルブは、その 2 つのポンプヘッドに含まれる 2 つだけであるため、障害発生の可能性も 抑えられます。ポンプのハードウェアに与える応力を削減するために、ピ ストンの加速と減速をゆっくりと行うことで衝撃を防止する円滑動作コン トロールを使用しています。これらの動きを制御できるように、高解像度 エンコーダユニットがポンプドライブに接続されています。このユニット は、回転を 65000 steps に設定し、各ステップのボリュームを約 300 pl に 変換します。



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の情報を示します。





設置について

機器が最適な性能で動作するためには、適切な環境に設置する必要があり ます。

電源について

モジュールの電源は、広範囲にわたる入力電圧に対応しており、『19ページ表1』に記載の範囲のいずれの入力電圧でも使用可能です。したがって、 モジュールの背面に電圧スイッチはありません。また、電源内に自動電子 ヒューズが装備されているため、ヒューズを外部に取り付ける必要はあり ません。

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。

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、カバーが開いていて、モジュール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。

警告
 モジュールの入力電圧が正しくありません
 装置を仕様よりも高い入力電圧に接続すると、感電の危険性や機器が損傷を受ける恐れがあります。
 → 使用するモジュールは、指定された入力電圧に接続してください。

注意

電源コネクタにが届くようにしてください。

緊急時に備えて、いつでも電源から装置を切り離せるようにしておく必要 があります。

- → 機器の電源コネクタは、簡単に手が届き取り外せるようにしておいてく ださい。
- → 機器の電源ソケットの後には、ケーブルを抜くために十分な空間を確保 してください。

電源コード

モジュールには、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。 どの電源コードの一方も、同じメス型です。電源コードのメス型側を、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コンセント合わせて設計 されています。

警告 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、感電や回路の短絡に至ることがあります。

- → 接地していない電源を使用して本装置を稼動しないでください。
- → また、使用する地域に合わせて設計された電源コード以外は、決して使用しないでください。

警告 指定外ケーブルの使用

アジレントが供給したものではないケーブルを使用すると、電子部 品の損傷や人体に危害を及ぼすことがあります。

→ 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。

警告提供された電源コードの目的外の使用

- 電源コードを目的外に使用すると、人体に危害を及ぼしたり、電子 機器に損傷を与えたりすることがあります。
- →この機器に付属の電源コードは、この機器以外には使用しないで ください。

設置スペース

モジュールの寸法と質量(『19 ページ 表 1』を参照)は、ほぼすべての机 やラボ作業台にモジュールを設置できるように設計されています。空気循 環と電気接続のために、本機器の両側に 2.5 cm (1.0 inches)、背面に 約 8 cm (3.1 inches)の空間が必要です。

作業台上に HPLC システム全体を設置する場合は、作業台がすべてのモジュールの質量に耐えるように設計されていることを確認してください。

モジュールは水平に設置して操作してください。

結露

注意

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	21.8 kg (48 lbs)	
寸法 (高さ × 幅 × 奥行き)	240 x 345 x 435 mm (9.3 x 13.5 x 17 インチ)	
入力電圧	100 - 240 VAC, ± 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	350 VA / 270 W / 922 BTU	最大値
使用周囲温度	4 – 55 ° C (41 – 131 ° F)	
保管周囲温度	-40 - 70 ° C ($-4 - 158$ ° F)	
湿度	$<$ 95 % r.h. \downarrow 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

タイプ	仕様	注釈
ハイドロリック システム	2 基のデュアルピストンを備 えた直列ポンプは弊社独自の サーボ制御可変ストロークデ ザイン。スムースな送液コン トロール。	
設定可能な流量 範囲	設定値 0.001 - 5 ml/min、 0.001 ml/min ステップ。	分解能 300 pl/step。
流量精度	≤0.07 % RSD または 0.005 min SD、いずれか大き い方 (0.2 -5.0 ml/min)。	一定の室温でのリテン ションタイムに基づい て。
流量真度	±1 % または 10 µ1/min、い ずれか大きい方。	脱気した H ₂ 0 を 10 MPa (100 bar) で送液。
圧力動作範囲	動作範囲最大 120 MPa(1200 bar)、最大 2 ml/min、5 ml/min で 80 MPa (800 bar)。	
圧力の脈動	<1 % または < 0.5 MPa (5 bar) の、いずれか大きい 方。	水 1 ml/min 送液時
圧縮率補正	溶媒選択による自動設定。	
グラジエント組 成	高圧バイナリ混合	

表 2 性能仕様

設置について 2 性能仕様

タイプ	仕様	注釈
ディレイボ リューム	Jet Weaver V35: <45 µl Jet Weaver V100: <75 µl	通常は JetWeaver を推 奨します。ディレイ ボ リュームが最小である必 要のあるアプリケーショ ンについては、流路から JetWeaver を取り外せま す。この場合、ディレイ ボリュームは、接続キャ ピラリのボリュームのみ で決定されます。
組成範囲	設定可能範囲: 0 - 100 %	推奨範囲:1 - 99 % ま たは、1 チャンネルにつ き 5 μ l/min、いずれか 大きい方。
組成精度	<0.15 % RSD、または 0.01 min SD、いずれか大きい 方。	0.2 -5.0 ml/min。一定 の室温でのリテンション タイムに基づいて
組成真度	± 0.35 % 絶対値(5 - 95 %、 0.2 - 5.0 ml/min)	水 / カフェイントレー サー
溶媒切り替えバ ルブ	デフォルト	ポンプの規準部品
一体型脱気装置	チャンネル数:2 チャンネルあたりの内部容 量:1.5 mL	
コントロール	Agilent LC ChemStation (B0.040.02 以降) EZChrom Elite (3.3.2 SP1 以 降) OpenLab (3.3.2 SP3) Masshunter (B.02.01 SP1 以 降)	

表 2 性能仕様

表 2 性能仕様

タイプ	仕様	注釈
ローカルコント ロール	Agilent インスタントパイ ロット (G4208A) (B.02.08 以降)	
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	
安全とメンテナ ンス	拡張診断機能、Agilent ラボ アドバイザによるエラー検出 と表示、リーク検出、安全 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 カシグナル。主要なメンテナ ンス領域における低電圧。	
GLP 機能	EMF (Early maintenance feedback)機能: 事前設定や ユーザー設定可能なリミット でのパージバルブ スイッチや ポンプの移動層のボリューム を継続的に追跡し、メッセー ジをフィードバック。メンテ ナンスとエラーの電子記録。	
ハウジング	全材料リサイクル可能。	



Agilent 1290 Infinity バイナリポンプ ユーザーマニュアル

。 モジュールの設置

3

モジュールの開梱 24
スタック構成の最適化 26
1 スタック構成 26
2 スタック構成 29
輸送用緩衝材の取り外し 31
ポンプの設置 32
ポンプへの配管 36
シールウォッシュオプションの取り付け 40

この章では、所有するシステムで推奨されるスタックのセットアッ プと、Agilent 1290 Infinity バイナリポンプの設置についての情 報を示します。



3 モジュールの設置 モジュールの開梱

モジュールの開梱

パッケージの不足および損傷

梱包箱の外観に破損などがある場合は、アジレントの営業所 / サービスオ フィスまで速やかにご連絡ください。サービス担当者に、機器が輸送中に 損傷を受けた可能性があることをご通知ください。

注意「到着時不良」の問題

モジュールに破損が見られる場合は、モジュールの設置を中止してください。機器の状態が良好であるか不良であるかを評価するには、アジレント による点検が必要です。

- → 損傷があった場合は、アジレントの営業およびサービスオフィスまでご 連絡ください。
- → アジレントのサービス担当者が、お客様の設置箇所における機器の点検 を行い、適切な初動動作を行います。

梱包チェックリスト

モジュール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。部品を識別するために、『「メン テナンス用部品と器材」191ページ図』の図解付き部品リストを確認して ください。不足または破損した部品があった場合は、Agilent Technologiesの営業およびサービスオフィスまでご連絡ください。

説明	数量
Agilent 1290 Infinity バイナリポンプユーザーマニュアル	1
Agilent 1290 Infinity システムマニュアル	1
Agilent Lab Advisor	1
LC HW ユーザー情報およびユーティリティ DVD	1
電源コード	1
溶媒キャビネットキット (ボトル 4 本付き)	1
アクセサリキット(『「アクセサリキット」209 ページ 図』を参照)	1
RRHD Eclipse Plus C18、2.1x50 mm,1.8 u	1

表 3 バイナリポンプ梱包明細

スタック構成の最適化

スタック構成の最適化

本モジュールを、Agilent 1290 Infinity 液体クロマトグラフの一部とし て使用する場合は、以下の構成で設置することで、最適な性能を得ること ができます。これらの構成でシステムの流路を最適化し、ディレイボ リュームを最小限に抑えます。

その他に可能なコンフィグレーションについては、『Agilent 1290 Infinity システムマニュアル』を参照してください。

1スタック構成

Agilent 1290 Infinity バイナリ LC システムの各モジュールを以下のコ ンフィグレーションで設置して、最適な性能が確実に得られるようにする 必要があります(『27ページ 図 2』と『28ページ 図 3』を参照)。このコ ンフィグレーションにより、ディレイボリュームを最小限に抑えるために 流路が最適化され、必要な設置スペースが最小になります。

Agilent 1290 Infinity バイナリポンプ バルブクラスタ コントロールは、 必ずスタックの最下部に設置する必要があります。

スタック構成の最適化



図 2 バイナリポンプを装備した 1290 Infinity の推奨スタックコン フィグレーション(前面図)

27

スタック構成の最適化



フィグレーション(背面図)

2 スタック構成

オート サンプラ用冷却モジュールがシステムに付いている場合は、重いモ ジュール(1290 Infinity ポンプとサーモスタット)を両方とも各スタッ クの最下部に設置して、スタックが高くならないようにした、2 スタック コンフィグレーションが推奨されます。オートサンプラ用冷却モジュール を追加しない場合でも、この構成を使ってスタックを低くすることが望ま しいことがあります。ポンプとオートサンプラ間には若干長いキャピラリ が必要になります(『29 ページ 図 4』および 『30 ページ 図 5』を参照)。



図 4 バイナリポンプを装備した 1290 Infinity の推奨 2 スタックコン フィグレーション(前面図)

スタック構成の最適化



AC 電源

図 5 バイナリポンプを装備した 1290 Infinity の推奨 2 スタックコン フィグレーション (背面図)

輸送用緩衝材の取り外し



3 モジュールの設置 ポンプの設置

ポンプの設置

- **必要な準備**: 設置スペースの決定 電源接続の準備 ポンプの開梱
 - 1 モジュールを水平な作業台の上に置きます。



2 モジュールの前面にある電源スイッチがオフになっている (スイッチが 飛び出ている) ことを確認します。

図 6 バイナリポンプの前面

- **3** Agilent 1290 Infinity バイナリポンプ背面の電源コネクタに電源ケー ブルを差し込みます。
- **4** 1290 Infinity バイナリポンプの背面に、必要なインターフェイスケー ブルを接続します。



電源コネクタ

図 7 バイナリポンプの背面

- 注記 Agilent 1290 Infinity システムでは、個々のモジュールを CAN ケーブルで接 続します。Agilent 1200 シリーズ インスタントパイロットは、どのモジュー ルの CAN バスにも接続できます。Agilent データシステムへの接続は、検出器 の内蔵 LAN ポートを通じて構築されます。全モジュールの中で最も取り込み速 度が速くなる検出器の LAN ポートを利用することをお勧めします。インスタン トパイロットまたは Agilent データシステムの接続についての詳細は、それ ぞれのユーザーマニュアルを参照してください。LAN アクセスの設定について は、『「LAN コンフィグレーション」251 ページ 図』を参照してください。
 - 5 モジュールの左下側にあるボタンを押して電源を ON にします。 電源ボタンは押し込まれた状態になり、ステータス LED が緑色になります。

注記 電源ボタンが飛び出た状態で、緑のランプが消えているときは、モジュールの 電源は切られています。

注記 モジュールは、デフォルトのコンフィグレーション設定で出荷されています。 これらの設定を変更するには、「8 ビットコンフィグレーションスイッチの設 定」のセクションを参照してください。



ポンプへの配管



ポンプは、溶媒切り替えバルブ、デガッサユニット、ポンプヘッド、圧力 センサ、パージバルブ、および Jet Weaver の間にチューブとキャピラリ 接続が設置された状態で出荷されます。この項では、追加の配管の設置に ついて説明します。
必要な部品:	番号	部品番号	説明
	1		他のモジュール
	1	G4220-68705	アクセサリキット
	1	G5067-1531	溶媒キャビネットキット

必要な準備: ポンプを LC システムに設置しておきます。

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。



1 前面カバーの両側にある止め具を押して、前面カバーを取り外します。

- **2** 1290 Infinity バイナリポンプが含まれるモジュールスタックの上に溶 媒キャビネットを置きます。
- 3 溶媒キャビネットの中にボトルを 4 本入れ、各ボトルにボトルヘッドア センブリをネジ留めします。
- 4 機器の左上隅にシャットオフバルブパネルを取り付けます。



 5 ボトルヘッドアセンブリから溶媒切り替えバルブのインレットコネクタ A1、A2、B1、B2 に溶媒チューブを接続します。水溶性溶媒(通常は チャンネル A1)には茶色のボトルを使用します。



ポンプへの配管

- 6 付属のシールを使用してチューブにラベルを貼り、溶媒キャビネットと 1290 Infinity バイナリポンプの止め具にチューブを固定します。
- **7** Jet Weaver の出口をオートサンプラに接続します。
- **8** 廃液チューブ(G4220-67000)を、ポート 5 および 6 のパージバルブの 出口に接続します。
- 9 波形廃液チューブをリークパネルの出口に接続します。
- 10 波形廃液チューブを廃液コンテナに挿入します。
- 11 ポンプ上部のモジュールのドレインチューブをポンプに通します。
- 12 初めて使用する場合には、システムをパージします(『「ポンプのプラ イミング」58ページ図』を参照)。

3 モジュールの設置

シールウォッシュオプションの取り付け

シールウォッシュオプションの取り付け



図 9 シールウォッシュ機能を使用したバイナリポンプ

1290 Infinity バイナリポンプには、オプションでシールウォッシュ機能 を使用できます。このオプションは、緩衝液を使用する場合や、ピストン やシールに堆積する可能性がある不揮発性溶媒または添加剤を使用する場 合にお勧めします。このオプションは、定期的にこれらの部品を自動でク リーニングするために使用します。

- **1** 洗浄溶媒ボトルを溶媒キャビネットに置きます。蒸留水とイソプロパ ノール (90/10) の混合液は多くのアプリケーションに適しています。
- 2 溶媒インレットチューブを溶媒ボトルに入れ、ボトルを閉じて、チューブをシールウォッシュポンプに接続します。
- 3 ウォッシュチューブの出口を廃液コンテナ中に挿入します。



Agilent 1290 Infinity バイナリポンプ ユーザーマニュアル

4 ポンプの使用

バイナリポンプの準備 42
HPLC システムでの藻の繁殖 43 藻の問題の予防 / 軽減方法 44
機器コントロール インターフェイスでのポンプの設定 45 概要 45 機器コンフィグレーション 46 ポンプユーザー インターフェイス (ダッシュボードパネ ル) 48 コントロール設定 51 メソッドパラメータの設定 53
ポンプのプライミング 58
溶媒情報 60

この章では、Agilent 1290 Infinity バイナリポンプの操作パラ メータについて説明します。



バイナリポンプの準備

バイナリポンプの準備

ポンプのベストパフォーマンスのために:

- 溶媒ボトルを入れた溶媒キャビネットは、必ずポンプの上(またはポンプより高い位置)に置いてください。
- 最適な性能のために、内蔵デガッサを使用します。0.5 mL/min より低い 流速とジェットウィーバー無しでのコンフィグレーションには必須で す。
- 真空脱気ユニット付きでポンプを使用するときは、ポンプを稼働する前に少なくともチャンネル当り5mLで脱気ユニットをフラッシュしてください。特にポンプシステムが一定時間(たとえば一晩中)電源が切られていた場合、そしてチャンネル内で揮発性溶媒の混合液が使用されている場合は、この作業が必要です。
- 溶媒フィルタが詰まらないように注意してください(溶媒フィルタなしでポンプを使用しないでください)。藻の繁殖を避けてください、 『「HPLC システムでの藻の繁殖」43ページ図』を参照。
- 定期的な間隔でポンプアウトレットフィルタとカラムフリットを点検し ます。詰まったポンプアウトレットフィルタは、その表面上の黒、黄、 緑様層で識別できます。
- 可能な限り溶媒チャンネルごとに最小流速 5 µL/min で送液し、未使用のポンプチャンネルにクロスフローしないようにしてください。
- バッファ溶液を使用した時は、システムの電源を切る前に水でフラッシュします。
- 濃度が 0.1 M 以上のバッファ溶液を長時間送液する時は、オプションの ウォッシュ用シール機能を使用してください。
- ・ 決して未使用のポンプをチャンネルに水を残したまま長時間(2~3
 日)放置しないでください。常に有機溶媒でフラッシュするか、または10%のイソプロパノールを水に加えてください。

ポンプの使用 4 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、誤った故障診断 やアプリケーション上のトラブルの原因となります。藻は水系溶媒中で繁 殖し、特に pH 4 ~ 8 の範囲で顕著です。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。

藻に関連する機器の問題

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。

- 溶媒フィルタの詰まりや、インレットあるいはアウトレットバルブへの 沈着が、不安定な送液、組成やグラジエントの問題、ポンプの故障を招 きます。
- インジェクタの前に装着する小さな孔径の高圧溶媒フィルタ等を塞ぎ、 システム圧が高くなります。
- PTFE フリットが微小成分により詰まり、システム圧の増大につながり ます。
- カラムフィルタを詰まらせ、システム圧を上昇させます。
- 検出器のフローセルウィンドウが汚れ、ノイズレベルが高くなります(検出器は流路の最後にあるモジュールのため、この問題が起こる頻度は 稀です)。

HPLC システムでの藻の繁殖

藻の問題の予防 / 軽減方法

- 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 水を使用してください。
- ・ 送液せず、移動相を数日間、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。
- ・ 茶色溶媒ボトル(溶媒ボトル、褐色(9301-1450)) 水系溶媒用、装置 に標準付属 - を使用します。
- 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 有機溶媒を水系溶媒に加えてください。

ポンプの使用 4 機器コントロール インターフェイスでのポンプの設定

機器コントロール インターフェイスでのポンプの設定

概要

以下のセクションで説明されるパラメータは、機器コントロール インター フェースで提供され、通常 Agilent 機器コントロール ソフトウェアを通 じてアクセスできます。詳細については、マニュアルやそれぞれのユー ザー インターフェイスのオンラインヘルプを参照してください。

機器コントロール インターフェイスでのポンプの設定

機器コンフィグレーション

機器コンフィグレーションダイアログボックスを使用して調べ、必要なら、 機器コンフィグレーションを変更してください。コンフィグレーション可 能なモジュールパネルは、コンフィグレーションに使用できる全てのモ ジュールのリストを含みます。選択したモジュールパネルは設定されてい るモジュールのリストを含みます。

自動コンフィグレーション:通信設定下では、ホスト名オプションまたは IP アドレス オプションを選択し、適切な値を入力して、ホストコン ピュータがハードウェア コンフィグレーションの自動検出できるようにし てください。システムが自動的に機器を設定し、それ以上のマニュアル コ ンフィグレーションは必要ありません。

バイナリポンプのコンフィグレーションパラメータは 2 つのセクションに あります:

- 通信
- ・ オプション

通信: このダイアログボックスでのパラメータは、オートコンフィグレー ションにより自動的に検出されます。

- デバイス名、
- ・タイプ ID、
- シリアル番号、
- ファームウェアリビジョン、
- ボタン 接続設定

機器コントロール インターフェイスでのポンプの設定

オプション:

・ 圧力単位:

ドロップダウンリストから圧力単位を選択してください (bar, psi また は MPa)。

・シールウォッシュ(オプション):

このチェックボックスをマークすると自動コンフィグレーション中にオプションのシールウォッシュを検出します。

• ISET (オプション):

このチェックボックスには、ISET が取り付けられていることを示す マークが付けられます。ISET コンフィグレーションをクリックして ISET Configuration ダイアログボックスを開きます。このダイアログ ボックスを使用して ISET エミュレーションのサンプラを設定できます。

溶媒タイプカタログを設定: 溶媒タイプカタログダイアログボックスを表示し、これにより溶媒キャリブレーションデータをインポートやエクス ポートできます。『「溶媒キャリブレーションテーブルのインポート」74 ページ 図』を参照してください。

詳細については、ユーザー インターフェイスのオンラインヘルプを参照し てください。

機器コントロール インターフェイスでのポンプの設定

ポンプユーザー インターフェイス (ダッシュボードパネル)

モジュールグラフィック

ポンプグラフィックの項目の意味と機能は以下のとおりです:

外部接点ボードがインストールされたことを示します。

ボトル中の溶媒レベルが緑のエリアで示されます。溶媒 レベルが指定の容量より少なくなるとそのエリアが黄に なり、ボトルが空になるとそのエリアは赤に変わります。 溶媒ボトルをクリックすると、溶媒ボトル容量 ダイアロ グボックスが表示されます。ボトルのツールチップは、 溶媒名を示します。

ISET オプションは据え付け済みだが、アクティブなメ ソッドがない (グレイ) か、インストール済みでアク ティブである (オレンジ) ことを示します (G4220A の み)。



溶媒切り替えバルブがあることを示します。グラフィッ クをクリックしてバルブを切り替えます。バルブが切り 替わるとアニメーションが表示されます。

100.00 bar

圧力設定値 赤線は現在の圧力上限を示し、緑のエリアは 現在の圧力を示します (テキストとしても示されます)。

現在の溶媒流速(mL/min)が、圧力表示上に表示されます。

機器コントロール インターフェイスでのポンプの設定

機器シグナル

以下のポンプシグナルが表示されます:

流量 現在の溶媒の流速 (mL/min)。

- **圧力** 現在のポンプ圧力 (bar、psi、MPa、『「機器コンフィグ レーション」46 ページ 図』を参照)。
- **圧カリミット** 現在の圧力上限
- 組成 A: B
 現在の溶媒組成を示します。溶媒切り替えバルブが取り 付けられると、チャンネルがグラフィックで表示されます。

機器コントロール インターフェイスでのポンプの設定

コンテキストメニュー

ダッシュボードパネルのコンテキストメニューは、以下のコマンドを含みます:

3	コントロール	ポンプの コントロール ダイアログボックス の表示。
Ś	メソッド	ポンプの メソッド設定 ダイアログボックス の表示。
	エラーメソッド を設定	現在使用しているメソッドでエラーが発生し た場合にロードされるメソッドを設定。
	装置の認識	モジュールの前部の LED が 2 ~ 3 秒点滅する 結果になります。
	ポンプのオン / オフ切り替え	ポンプのオン / オフのステータス切り替え。
	溶媒ボトル容量	ポンプの 溶媒ボトル容量 ダイアログボック スの表示。
	パージオン / オ フ	システムのパージをコントロール。
	プライミング オ ン / オフ	最初にポンプヘッドに溶媒を引き込むための プライミングを実施。
	コンディショニ ング オン / オフ	ポンプコンディショニングのオン / オフ切り 替え。コンディショニング機能は、ポンプの 流路内の気泡を除去するのに有効。

機器コントロール インターフェイスでのポンプの設定

コントロール設定

バイナリポンプのコントロール パラメータは6つのセクションに分かれま す:

- ・ ポンプ
- ウォッシュ用シール
- ・ 自動オン
- ・パージ
- ・ プライミング
- ・ コンディショニング
- 表 4 ポンプコントロール パラメータ

パラメータ	範囲	説明
ポンプ		ポンプのスイッチを オン、オフ または スタンバイ の状態にできます。 スタンバイ の状態では、ポンプのモーターはまだアク ティブで、ポンプのスイッチを再び入れるとき再度初期 化する必要はありません。
ウォッシュ用 シール		 ウォッシュ用シールを一度実行(シングル洗浄)または 周期的(周期的)に設定できます。 オフ:ウォッシュ用シールを使用しません。 シングル洗浄:シールは指定時間パージされます。 周期的:設定された時間(分)周期的洗浄を適用します。 このオプションは、ポンプがシールウォッシュ機能を有する場合のみ使用できます。シールウォッシュ機能はモジュールが認識しています。もしウォッシュ用シールがインストールされている場合は、主要なシールの寿命を長くするため使用することが推奨されます。
ウォッシュ用 シール ラン モード		 ウォッシュ用シールを使用する時を設定できます。 オフ:ウォッシュ用シールは非アクティブです。 ポンプオンの時、オン:ウォッシュ用シールは、ポン プがオンの時のみアクティブです。 常時オン:ウォッシュ用シールは、ポンプがオンまた はスタンバイ モードの時にアクティブです。

機器コントロール インターフェイスでのポンプの設定

表 4 ポンプコントロール パラメータ

パラメータ	範囲	説明
自動オン		指定した日付 / 時間にモジュールをオンにすることがで きます。この機能は、モジュール電源スイッチがオンの 場合のみ使用できます。
パージ	時間:0- 100.00 min0.01 ステップ。 流量:各チャン ネルについて 0.000 - 5.000 mL/min、 0.001 ステップ (最大 10.000 mL/min)。	 パージ パラメータの設定とアクティベーション 自動パージバルブは、システムをパージするために使用できます。 そのプロセスは、使い易くするため自動化されました。 オフ:パージをオフにします。 オン:装置はパージされます。 パージ中のパージ流量、時間、組成を設定する必要があります。 パージ時間が終了するとすぐに、モジュールは自動的に分析コンディションに再度切り替わります。
プライミング		プライミングを開始するにはオンを、オフにするにはオ フを選択します。 プライミング機能は、空の溶媒ラインを充填したり、空 気がポンプヘッドに侵入した場合に役立ちます。モ ジュールは 4 つのポンプドライブを同時に使って高速で 溶媒を吸引し、それを自動パージバルブの廃液ポジショ ンに供給します。これを 20 回実行して、プロセスは終了 します。
コンディショニ ング	少なくとも 200 bar (> 500 bar がよ り良い)。	もし超過圧力または組成リップルを検出した場合はこの 機能を使用して、溶媒タイプ(水/有機溶媒、または特 殊溶媒/溶媒混合)が正しくセットされており、ポンプ に漏れの徴候がないことを確認してください。 例えば、溶媒が切れた後、長いスタンドバイ時間の後、 またはサービスあるいは修理の後で、もしポンプが空気 を含んでいる場合、コンディショニングが必要になるこ とがあります。

機器コントロール インターフェイスでのポンプの設定

メソッドパラメータの設定

バイナリポンプのメソッド設定パラメータは、8つのセクションに分かれます:

- 流量
- 溶媒 A と B
- ・ ストップタイム
- ・ ポストタイム
- 圧力リミット
- ・ タイムテーブル
- 詳細
- 外部接点
- 表 5 メソッド パラメータ

パラメータ	範囲	説明
流量	0.00 - 5.00 mL/min、 0.001 ステッ プ。 推奨流量範囲: 0.05 - 5.00 mL/min.	流量は、カラムに沿って溶離液が移動する速度です。正確なリテンション タイムとピーク測定を確実にするためには、流量を一定に保つことが重要です。ポンプ システムが部分的失敗、ポンプ システム中の空気、移動相の粘性変化、または温度変化の結果、流量の変動が生じます。
溶媒 A と B		各チャンネルに対して、2 種類の溶媒のどちらを供給す るかを選択できます。溶媒 B のパーセンテージは、0 から 100 % までの任意の値に設定できます。溶媒 A は、 常に残りの容量を送液します (100 - %B)。溶媒 B の チェック ボックスで、溶媒チャンネルをオン (チェック あり)またはオフ (チェックなし)にできます。 圧縮率 セクションの 溶媒タイプの使用 チェックボック スがオンのとき (『「詳細設定」55 ページ 図』を参照)、 下向き矢印をクリックして 一般的 溶媒またはキャリブ レーション済みの 溶媒 を選択します。

機器コントロール インターフェイスでのポンプの設定

表 5 メソッド パラメータ

パラメータ	範囲	説明
ストップタイム	0.01 - 99999 min または インジェ クタと同様 / 無 制限 (分析時間 無限)。	タイムリミットが、分析用の時間制限を設定します。ス トップタイムの後、全てのグラジエントが停止し、ポン プ パラメータは初期値に戻ります。ポンプは、完全な分 析システムのストップタイムマスターとして使用できま す。検出器のストップタイムが 無制限 に設定さ r てい てもポンプが検出器を停止させます。無制限の場合、メ ソッドをマニュアルで停止する必要があります。
ポストタイム	0.01 - 99999 min また は オフ (0.0 min)。	機器は、次の分析開始を遅らせるため、ポストタイム中 はノットレディの状態のままです。ポストタイム を使用 して、溶媒組成変更後カラムが平衡化できるようにしま す(例えば、グラジエント溶出後)。
圧カリミット	上限: 120 MPa (1200 bar、 17400 psi)最 大流速2mL/min まで。2mL/min から5mL/min で、80 MPa (800 bar、 11600 psi)まで 一定比率で。 下限:0以上、 圧力リミット上 限の間の任意の 数値	 ポンプの最大圧カリミットと最小圧カリミットを設定してください。 上限は、ポンプがそれ自体でオフに切り替わり、超過圧力に対して分析システムを保護する圧力上限です。 下限は、例えば溶媒ボトルが空の場合、ポンプ自体でオフに切り替わる圧力下限であり、空気が送られてシステムが損傷してしまうことを避けます。
タイムテーブル		を参照 『「タイムテーブル設定」57 ページ 図』
詳細		『「詳細設定」55 ページ 図』を参照
外部接点		 外部接点セクションにより、外部接点の切り替えを設定できます。 注記 外部接点セクションは、BCD/外部接点ボードが据付済みの場合のみあります。

機器コントロール インターフェイスでのポンプの設定

詳細設定

バイナリポンプの詳細メソッド設定パラメータは、3 つのセクションにあります:

- ・ 最小ストローク
- 丘縮率
- ・ 最大流量 グラジエント

表 6 詳細メソッド パラメータ

パラメータ	範囲	説明
最小ストローク	20 - 100 µL	ストローク容量は、モジュールのパフォーマンスとシー ルの寿命間を最適化するために使用します。パフォーマ ンスのためには、低ストローク容量の方が、脈流を低く 抑えることが可能ですが、ストローク容量を大きくする ことでポンプシールの寿命を伸ばすことができます。 もし自動がアクティブの場合、、ポンプはジェットウィ バーの位置を考慮して最適なストロールボリュームを設 定します。 同期: 両方のチャンネルのストロークを同期させるには、 このオプションを選択します。チャンネル B の値は、 チャンネル A の値と同じに設定されます。こうすること によってフローティング障害が機器の性能に影響しない ようにします。

機器コントロール インターフェイスでのポンプの設定

表 6 詳細メソッド パラメータ

パラメータ	範囲	説明
圧縮率		 移動相の圧縮率は、ポンプのパフォーマンスに影響します。最適な流量真度と混合パフォーマンスを得るために、使用されている移動相に従ってパラメータを設定できます。 溶媒タイプの使用: このチェックボックスをオフにすると、圧縮率フィールドが表示され、圧縮率を入力できます。 チェックボックスが選択(オン)されていると、圧縮率フィールドは非表示になり、拡張圧縮率キャリブレーションが有効になります。溶媒 セクション内のコンボボックスを使用して、ドロップダウンリストから必要なキャリブレーション済みの溶媒を選択してください。
最大流量 グラジ エント	1.000 - 1000.000 mL/min /min ステップ 0.001 mL/min/mi n 初期値: 100.000 mL/min/ min	溶媒フローの変更速度にリミットを設定して、分析カラ ムを保護できます。 流量上昇と流量下降 用に、個々の値 を設定できます。

機器コントロール インターフェイスでのポンプの設定

タイムテーブル設定

タイムテーブルの、**タイム**フィールドに時間を入力し、続くタイムテー ブルのフィールドに適切な値を入力して、分析中ポンプパラメータの変更 をプログラムしてください。流速は、ゼロ時間または最後に変更を設定し た時間から直線的に変化します。他のパラメータは、タイムテーブルで設 定された時間に即座に変わります。

詳細タイムテーブル表示が、タイムテーブル表示を標準モードと詳細モー ド間で切り替えます。

以下のパラメータが変更可能です:

- 接点の変更
- 流量変更
- ・ 上限圧カリミット変更
- 溶媒組成変更 溶媒セクションで有効化した溶媒のみを使用できます。
- 機能中心表示 このチェックボックスにより、タイムテーブルの代わり にパラメータ変更内容を表示できます。

ポンプのプライミング

ポンプのプライミング

溶媒が交換されたり、また一定時間(たとえば一晩中)ポンプシステム の電源を切った場合、溶媒リゾルバ、真空脱気ユニット(システムで使用 できるとき)、ポンプ間の溶媒チャンネル中に気泡が再度拡散します。揮発 性化合物を含む溶媒は、化合物が揮発し、濃度が変化します。そのため、 アプリケーションを開始する前に、ポンプシステムのプライミングが必要 です。

- **1** パージ流量をチャンネル当り 3 5 ml/min に設定して、制御ソフト ウェアのパージを開始してください。
- 2 少なくとも 30 ml の溶媒を使用して、すべてのチューブを洗浄します。

表 7 さまざまな目的に対するプライミング用溶媒の選択

内容	溶媒	注釈
インストール後	イソプロピルア ルコール	システムから気泡を洗い出すために最 適な溶媒
逆相と順相を切り替える際(両方と も)	イソプロピルア ルコール	イソプロピルアルコールは、順相溶媒 と逆相溶媒の両方に混和性がありま す。
インストール後	エタノールまた はメタノール	イソプロピルアルコールが入手できな い場合の代用 (第 2 の選択肢)
緩衝液使用中にシステムを洗浄する	再蒸留水	析出した緩衝液の再溶解に最適な溶媒
溶媒を交換した後	再蒸留水	析出した緩衝液の再溶解に最適な溶媒
システムを長時間オフにする前に	有機溶媒または 10 % イソプロピ ルアルコール水	

注記

空のチューブの呼び水はポンプを使用して行わないでください(ポンプは、乾燥させないでください)。ポンプを使った呼び水を行う前に、シリンジを使用して溶媒を十分に吸引し、ポンプインレットへのチューブを完全に溶媒で満たしてください。

ポンプのプライミング

もしシステムをドライで運転したり、または空気がポンプ内に拡散した 場合、空気を再度除去するための追加手順が必要になることがありま す。以下の手順に従うとベストで最速の結果が得られます。

- 溶媒を両方のチャンネルについてイソプロピルアルコールに変更してく ださい。
- 2 プライム機能をオンにします。
- **3** システムを、10 ml(混合比 50/50) で 10 min.
- 4 イソプロピルアルコールに適したカラムを接続し、最大圧力限界値をカ ラムの限界値に設定します。
- 5 混合比 50/50、カラムの限界値に近い圧力を生成する流量でシステムを 実行します。
- 6 圧力の変動を観察します。圧力が安定すると、システムは脱気されてい ます。
- 7 分析条件に従って溶媒とカラムを変更し、システムをパージして溶媒を 交換します。

4 ポンプの使用 溶媒情報

溶媒情報

概要

溶媒を使用するときは、次の注意に従ってください。

- 藻の増殖を避けるための推奨事項に従ってください。『「HPLC システム での藻の繁殖」43ページ 図』を参照してください。
- 小さな粒子がキャピラリとバルブを永久的にブロックすることがあります。従って、常に 0.4 µm フィルタで溶媒を必ず濾過してください。
- 流路内の部品の腐食の原因となる溶媒の使用は避けるか、最小限にしてください。フローセルやバルブ材などの異なる材料に対して示されたpH範囲に関する仕様や、以降のセクションの推奨事項を考慮してください。

流路内の材質

以下の材質がこのモジュールで使用されています:

部分	材質
デガッサチャンバ	TFE/PDD 共重合体、FEP、PEEK、PPS
マイクロフルイ ディック構造 ¹	SST
SSV	PEEK, FEP, PFA, A1 ₂ 0 ₃ -ベースのセラミック製、ル ビー、サファイア、SST
パッシブインレッ トバルブ	SST、金、ルビー、Zr0 ₂ -ベースのセラミック製、タン タル
アウトレットバル ブ	SST、金、ルビー、ZrO ₂ ベースのセラミック、タンタ ル
ポンプヘッド	SST
ピストン	ZrO ₂ ベースのセラミック
ピストン / ウォッ シュ用シール	UHMW-PE, SST
圧力センサー	SST
自動パージバルブ	ポリイミド、SST、DLC

¹ ジェットウィーバー、熱交換器



材質情報

流路内の材質は、数十年かけて HPLC 分析用に高品質の設備を開発してきた Agilent 社の経験に基づき、注意深く選択されています。これらの材質は、 標準の HPLC 条件下で優れた堅牢性を発揮します。特殊な条件については、 材質情報セクションを参照するかまたは Agilent 社に連絡してください。

免責条項

次のデータは外部の情報源から収集したもので、参考としてお使いいただ くためのものです。Agilent 社は、このような情報の完全性や正確性を保 証することはできません。データは、適合性ライブラリに基づいており、 それは UHPLC システム、溶媒、溶媒混合液、サンプルの、特殊であっても 非常に多様な条件下での長期間の寿命を推定するために固有のものではあ りません。情報は、金属イオン、錯化剤、酸素などのような不純物の触媒 効果により、一般化できません。純粋化合物の腐食以外に、電食、静電気 的変化(特に非伝導性有機溶媒)、ポリマの膨張など他の影響も考慮する必 要があります。得られたデータの大部分は室温(通常 20 - 25 °C、68 -77 °F)に言及しています。腐食の可能性がある場合、温度が上昇すると 一般に腐食が加速します。疑わしいときは、材料の化学的適合性について の技術文献を参照してください。

PEEK

PEEK (ポリエーテル-エーテル ケトン) は、耐薬品性、機械的安定性、 熱安定性に関して優れた特性を併せ持っています。 pH1 ~ 12.5 の範囲 で安定であり、一般的溶媒の多くに対し不活性です。次のような化合物と の不適合性についてもいくらか知られています。クロロホルム、塩化メチ レン、THF、DMSO > 1 %、強酸(硝酸 > 10 %、硫酸 > 10 %、トリクロロ 酢酸、スルホン酸)、ハロゲン、または含水ハロゲン溶液、フェノール、誘 導体 (クレゾール、サリチル酸など)。室温を超える状態で使用すると、 PEEK は塩基やさまざまな有機溶媒に反応し、膨張を引き起こすことがあり ます。

ポリイミド

Agilent 社は、バルブ内のロータシールとオート サンプラのニードルシートに、半結晶性ポリイミドを使用しています。ポリイミドの供給元のひとつとして DuPont 社があり、そのブランド名は Vespel で、Agilent 社でも使用されています。

ポリイミドは、ほとんどの有機溶媒で、pH1 ~ 10 の 範囲で安定です。濃 縮鉱酸(例、硫酸)、氷酢酸、DMS0、THF とは適合しません。アンモニア (例、塩基性条件下でのアンモニウム塩)のような求核性物質または酢酸に よって分解します。

ポリエチレン(PE)

Agilent 社は、1290 Infinity ポンプで使用される黄色のピストンと ウォッシュ用シール用に、そして 1260 Infinity ポンプの順相アプリケー ション用に、UHMW (超高分子量)-PE/PTFE 混合物を使用しています。

ポリエチレンは、pH1 ~ 12.5 の 範囲の酸と塩基を含め、大抵の一般的無 機溶媒に対し、高い安定性をもちます。メタノール、アセトニトリル、イ ソプロパノールなど、クロマトグラフシステムで使用される多くの有機溶 媒に適合します。脂肪族、芳香族、ハロゲン化炭化水素、THF、フェノール や誘導体、濃酸と濃塩基とでは、安定性が限定されます。順相アプリケー ションに対しては、最大圧力は 200 bar に限定されます。

タンタル(Ta)

タンタルは、大抵の一般的 HPLC 溶媒と、無水フッ酸と三酸化硫黄をもつ 酸を除くほぼ全ての酸に対して不活性です。強塩基(例、水酸化物溶液 > 10 %、ジエチルアミン)により腐食します。無水フッ酸やフッ化物との使 用は推奨されません。

ステンレススチール (SST)

ステンレススチールは多くの一般的な溶媒に対して不活性です。1 から 12.5 の pH 範囲の酸や塩基の存在下では安定です。pH 2.3 より低い酸に より腐食することがあります。以下の溶媒によっても腐食することがあり ます。

- ハロゲン化アルカリ化合物およびその酸溶液(例、ヨウ化リチウム、塩 化カリウムなど)と、ハロゲン化物の水溶液。
- ・ 硝酸、硫酸などの高濃度の無機酸、特に高温の有機溶媒(クロマトグラフィーメソッド上可能であれば、ステンレスに対する腐食性の低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。

 $2 \text{ CHC1}_3 + 0_2 \rightarrow 2 \text{ COC1}_2 + 2 \text{ HC1}$

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。

- ・ 過酸化物(THF、ジオキサン、ジイソプロピルエーテルなど)を含む可能
 性がある、クロマトグラフクラスのエーテル。このようなエーテルは、
 過酸化物を吸着する乾性アルミニウム酸化物を使用して濾過してくださ
 い。
- 有機溶媒中の有機酸溶液(酢酸、ギ酸など)。例えば、酢酸1%のメタノール溶液は鋼鉄を腐食します。
- ・ 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。

ダイアモンド様炭素 (DLC)

ダイアモンド様炭素は、ほぼ全ての一般的な酸、塩基および溶媒に対して 不活性です。HPLC アプリケーションに対する不適合は報告されていませ ん。

フューズドシリカとクォーツ (SiO₂)

フューズドシリカは、1290 Infinity フロー セルとキャピラリに使用され ています。クォーツは、旧モデルのフローセル窓用に使用されています。 フッ化水素酸とフッ化物を含む酸性溶媒を除く全ての溶媒に対し、不活性 です。強塩基によって腐食するため、室温で pH 12 を超える状態で使用し ないでください。フローセル窓が腐食すると、測定結果に悪影響を及ぼす ことがあります。pH が 12 を超える場合、サファイアウィンドウ付きのフ ロー セルを使用することを推奨します。

金

金は指定の pH 範囲内では、一般的な全ての HPLC 溶媒、酸および塩基に 対して不活性です。錯化性シアン化物や王水のような濃酸により腐食する ことがあります。

ジルコニア 酸化物 (ZrO₂)

ジルコニア酸化物は、ほぼ全ての一般的な酸、塩基および溶媒に対して不 活性です。HPLC アプリケーションに対する不適合は報告されていません。

プラチナ / イリジウム

プラチナ / イリジウムは、ほとんどすべての一般的な酸、塩基および溶媒 に対して不活性です。HPLC アプリケーションに対する不適合は報告されて いません。

フッ素化ポリマ(PTFE、PFA、FEP、FFKM)

PTFE (ポリテトラフルオロエチレン)、PFA (パーフルオロアルコキシ)、 FEP (フッ素化エチレン プロピレン)などのフッ素化ポリマは、ほぼ全て の一般的な酸、塩基、溶媒に対し不活性です。HPLC アプリケーションに対 する不適合は報告されていません。FFKM は過フッ素化ラバーであり、ほと んどの化学物質に対しても耐性を示します。エラストマであることから、 ハロゲン化炭化水素などの有機溶媒中で膨張することがあります。

サファイア、ルビー、Al₂03-ベースのセラミック製

サファイア、ルビー、アルミニウム酸化物 Al₂0₃ ベースのセラミック製 は、ほぼ全ての一般的な酸、塩基、溶媒に対し不活性です。HPLC アプリ ケーションに対する不適合は報告されていません。





Agilent 1290 Infinity バイナリポンプ ユーザーマニュアル

性能の最適化

5

ディレイボリュームとエクストラカラムボリューム 68 ディレイボリューム 68 最適ディレイボリュームの設定方法 69 高分解能を達成する方法 71 溶媒キャリブレーションテーブルの使用 74

この章では、追加デバイスの性能または使用の最適化のヒントについて説明します。



5 性能の最適化

ディレイボリュームとエクストラカラムボリューム

ディレイボリュームとエクストラカラムボリューム

ディレイボリュームは、ポンプ内のミキシングポイントとカラム入り口間の システムボリュームと定義されます。

カラム外ボリュームは、カラム内のボリュームを除外した、注入ポイントと 検出ポイント間のボリュームと定義されます。

ディレイボリューム

グラジエント分離では、このボリュームによって、ポンプ内で変化してい る混合物間でディレイが生じるため、カラムへの到達が変化します。ディ レイは、流速とシステムのディレイボリュームに依存します。これは、各 HPLC システム内では、各分析の開始時に、グラジエントプロファイル内に 追加アイソクラティック溶出の時間が存在することを意味します。通常、 グラジエントプロファイルは、ポンプでのグラジエントプログラムを使っ てレポートされますが、ディレイボリュームは、クロマトグラフィに影響 するにもかかわらず、レポートには含まれません。この効果は、低流速と 小さなカラムボリュームでは影響が大きくなり、グラジエントメソッドを 移管した場合に、、最大の影響を与えます。そのため、高速グラジェント分 離の場合、小さなディレイボリュームにすることが重要です。特に、質量 分析検出にしばしば使用されるナロー ボアカラム (例: 2.1 mm 内径)の 場合は、注意する必要があります。

システム内のディレイボリュームには、ポンプ内のミキシングポイントからのボリューム、ポンプとオートサンプラ間の接続、オートサンプラを経由する流路のボリューム、オートサンプラとカラム間の接続が含まれます。

最適ディレイボリュームの設定方法

ポンプの物理的なディレイボリュームは、主として、Jet Weaver ミキサー の使用に依存します。UV 検出の場合は Jet Weaver を使うことが必須です が、質量分析検出の場合はディレイボリュームを削減するために Jet Weaver を使わないことを選択できます。これが意味を持つのは、超高速グ ラジエント操作(0.5 min以下)または非常に小さなボリュームのカラム を使う場合だけです。Jet Weaver を迂回した場合は、オートサンプラへの 接続チューブが、パージバルブから直結されることになります。

注記

流路から Jet Weaver を切断する前に、有機溶媒でフラッシュします。水や緩 衝液が Jet Weaver 内部に残らないようにしてください。残っていると藻やバ クテリアなどの微生物が繁殖する場合があります。

ポンプ内のディレイボリュームを増やすことが必要な場合があります。具体的に言えば、UV 検出を使う場合に、UV の吸収率が非常に高い化合物を 移動相に追加する場合です。この場合は、どんなポンプノイズでも拡大さ れます。最も一般的な例は、蛋白質とペプチドの分析でトリフルオロ酢酸 (TFA)を使うことです。この影響は、ミキサー容量を増加させることで緩 和できます。

以下のような異なる Jet Weaver の設定が利用可能です。

品目 Jet Weaver 35 µL/ 100 µL (G4220-60006) には、同一ユニット内に2つの代替ボリュームがあります。

低ボリューム 35 μ 1 から高ボリューム 100 μ 1 への切り換えは、アンイ ンストール、前面から背面への回転、そして再インストールすることに よって行います。『「構成の変更または Jet Weaver の交換」146 ページ 図』を参照してください。ミキシングボリューム (結果としてはディレ イボリューム) は、65 μ 1 だけ増加し、TFA のような添加剤を使った場 合のベースラインパフォーマンスが改善されます。Jet Weaver の構成 は、添付されている RFID タグによって自動的にログに記録されます。 5 性能の最適化 最適ディレイボリュームの設定方法

380 µL Jet Weaver 高性能ミキサは、例えば、トリフルオロ酢酸(TFA)のような高い吸光度を持つ移動相を使用したり、A チャンネルと B チャンネルで UV 吸収の異なる移動相を混合するといったアプリケーションに特化して使用をお勧めします。

ポンプで作成された溶媒パッケージは、溶媒が検出器フローセルに達す るまで持続します。吸収変動はベースライン ノイズとして現れ、混合ノ イズとも呼ばれます。不純物質定量または最低レベルの化合物検出のよ うなアプリケーションでは、ノイズを最小に抑える必要があります。 380 µL Jet Weaver は、混合をかなり改善するので、ベースライン ノイ ズを減らし、検出感度を向上させます。特許の Agilent マイクロフルイ ディック技術によって、380 µL という低内部ボリュームで高い混合パ フォーマンスが得られます。

高分解能を達成する方法

分離分解能を向上させると、定性的および定量的分析機能が改善され、分離可能なピークが増え、分離をさらに加速する可能性が増えます。このセクションでは、以下の要素を検討することによって、分離度を向上させる 方法について説明します。

- 選択性の最適化
- 小さな粒子径の充填剤
- 長いカラム
- 浅いグラジエント、高流量

2 つのピーク間の分解能は、次の分解能方程式で表現できます。

$$Rs = \frac{1}{4}\sqrt{N}\frac{(\alpha - 1)}{\alpha}\frac{(k_2 + 1)}{k_2}$$

変数の意味は次のとおりです。

- R_s= 分離度、
- N = 理論段数(カラム効率の目安)、
- a= 選択性(2 つのピーク間)、
- k₂=2 番目のピークの保持係数(旧称、キャパシティ係数)

分離度に大きく影響する項は選択性、a であり、この項を変更する要素に は、固定相のタイプ (C18、C8、フェニル基、ニトリルなど)、移動相、そ して分離対象の溶質間の選択性の差異を最大化するための温度などの変更 が含まれます。この操作は分析の本質的な部分であり、最も効率的に行う には、自動メソッド開発システムを使います。このシステムを使えば、異 なるカラムや移動相の広範囲の条件を順序付けられたスカウティングプロ トコルにより評価できます。このセクションでは、より高い分離度が得ら れる固定相や移動相の選択について説明します。自動メソッド開発システ ムを移動相選択のために利用し、スカウティングの各ステップで短いカラ ムを利用することで、高速分析を行うことができます。

分離能方程式によれば、次に重要な項は理論段数、言い換えると、効率 N であることがわかります。この項はいくつかの方法で最適化できます。N

は粒子径に反比例し、カラムの長さに直接比例します。したがって、小さ な粒子径と長いカラムを選択すれば、段数が大きくなります。圧力は、粒 子径の逆二乗で増加し、カラムの長さに比例します。これが 1290 Infinity LC システムが 120 MPa (1200 bar) までの圧力とサブ 2 µm の 粒子に対応し、カラムの長さを 100 mm または 150 mm まで伸ばせるよう に設計された理由です。また、100 mm と 150 mm のカラムを接続して 250 mm の長さに構成することもできます。分離能は N の平方根で増加す るため、カラムの長さを倍にすれば、分離度は 1.4 倍増加することになり ます。達成可能な数値は、移動相の粘性に依存します。これは圧力に直接 関係するためです。メタノール混合液は、アセトニトリル混合液よりも高 い背圧を生成します。アセトニトリルは粘性が低いことに加えて、ピーク の形状が優れていて狭いため、高い頻度で使用されますが、選択性はメタ ノールの方が優れています(特に、500 Daより小さい分子の場合)。粘性 は温度を上げれば減らせますが、その場合は分離の選択性が変化する可能 性があることに注意してください。経験によって、温度を上げたり下げた りすることでどのように選択性が変化するかが分かってきます。流量や圧 力を上げるとカラム内の摩擦熱が増加するため、拡散が増加し、選択性を 若干変化させる原因となることに注意してください。この増加や変化は、 分離度が減少することで確認できます。後者の場合は、サーモスタットの 温度を数度下げれば相殺できる可能性があります。これも経験で正解を導 くことができます。

Van Deemter 曲線は、STM カラム内の最適流速が、大きな粒子に対しては 多く、流速が増大するにつれて平準化していくことを示しています。STM カラムの最適に近い標準流速は:内径 4.6 mm カラムでは、2 ml/min 内径 2.1 mm カラムでは 0.4 ml/min です。

イソクラティック分離では、保持係数 k を増加させると、溶質の保持時間 が増加するため分離度が向上します。グラジエント分離では、リテンショ ンは次の式の k* で表現されます。

$$k^* = \frac{t_G}{\Delta\%B} \cdot \frac{F}{V_m} \cdot \frac{100}{S}$$

変数の意味は次のとおりです。

- k* = 平均 k 值
- ・ $t_G = グラジエントの時間の長さ(または、グラジエントのセグメント)$ (分)
- F = 流量(m1/分)
- V_m = カラムディレイボリューム
- D%B = グラジエント中の溶媒 B が変化する割合
- S = 定数 (分子が小さい場合は、約 4 ~ 5)

この方程式によれば、k、ひいては分離度は、グラジエントを浅く(目安としては、2から5%/minの変化)、流速を上げ、ボリュームカラムを小さくすれば、増加することがわかります。流量を倍にしてもグラジエント時間を半分にすれば、k^{*}を変化させずに分離性能を維持できて、時間は半分で済むようになります。最近発表された研究には、短いSTMカラムを(40°Cより上の温度で)高速で動作させることによって、長いSTMカラムよりも高いピークキャパシティを生成する方法が示されています。(Petersson et al., J. Sep. Sci, 31, 2346-2357, 2008, Maximizing peak capacity and separation speed in liquid chromatographyを参照)。

5 性能の最適化

溶媒キャリブレーションテーブルの使用

溶媒キャリブレーションテーブルの使用

溶媒キャリブレーションテーブルのインポート

RC.NET ベースの Agilent グラフィック ユーザー インターフェイス (ChemStation、EZChrom Elite、OpenLab など)は、HPLC でごく一般的に 使用される溶媒についてのデータが組み込まれています。このデータには 溶媒特性が含まれており、最適な流量を組成真度を得るためのコポンプコ ントロールの最適化に使用されます。

ソフトウェアに、使用する溶媒が含まれていない場合は、Agilent 社の ウェブサイト http://www.chem.agilent.com/en-US/Support/Downloads/firmware/Pages/LC.aspx で、追加ライブラリをチェックしてくださ い(登録が必要です)。そのライブラリでは、更新された最適なデータが提 供されています。

使用する溶媒がユーザー インターフェイスでもライブラリでも見つからない場合は、一般的な溶媒を使用してください。100%から 50 % までの水溶液は純水と特性が同等であるため、"一般的な溶媒"を選択することで、良好な結果を得ることができます。有機溶媒の割合が高い溶媒の場合は、"一般的な有機溶媒"を選択してください。

性能の最適化 5

溶媒キャリブレーションテーブルの使用

ChemStation での溶媒キャリブレーションのインポート

- **1** メニュー機器 > 機器コンフィグレーションに移行します。
- 2 機器コンフィグレーション スクリーンで、使用するモジュールを選択 し、設定をクリックします。
- 3 溶媒タイプカタログを設定をクリックしてください。
- **4 溶媒タイプカタログ** で、**インポート**をクリックします。
- 5 溶媒キャリブレーション テーブルの場所にナビゲートし、開くをク リックします。
- 6 ここで、新しい溶媒が溶媒タイプカタログに現れます。
- これでインポートされた溶媒を、溶媒タイプの選択に使用できます。
 『53ページ表5』を参照してください。

5 性能の最適化

溶媒キャリブレーションテーブルの使用

Agilent 1290 Infinity バイナリポンプ ユーザーマニュアル



6 トラブルシューティングおよび診 断

モジュールのインジケータとテスト機能の概要 78
ステータスインジケータ 79
電源インジケータ 79
モジュールステータスインジケータ 80
使用可能なテストとユーザーインタフェース 81
Agilent Lab Advisor ソフトウェア 82

トラブルシューティングと診断機能についての概要



モジュールのインジケータとテスト機能の概要

モジュールのインジケータとテスト機能の概要

ステータスインジケータ

モジュールには、モジュールの稼動ステータス(プレラン、ラン、エラー 状態)を示す 2 つのステータスインジケーターが装備されています。ス テータスインジケーターによって、モジュールの動作状態を一目で確認す ることができます。

エラーメッセージ

モジュールの電子、機械、または流路系統に障害が発生した場合は、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。各メッセージ について、障害の簡単な説明、その原因、および対策を示します(「エ ラー情報」の章を参照)。

テスト機能

トラブルシューティングと内部部品交換後の動作確認のために、一連のテ スト機能が用意されています(「テスト機能とキャリブレーション」を参 照)。

診断シグナル

モジュールには、問題の診断に使用できるシグナルがいくつかあります(ランプの内部温度、電圧、電流)。これらは、Agilent ChemStation ソフ トウェアに標準シグナルと同様に追加できます。

トラブルシューティングおよび診断 6 ステータスインジケータ

ステータスインジケータ

モジュールの前面には、2 つのステータスインジケータがあります。左下 のインジケータは電源ステータスを示し、右上のインジケータはモジュー ルステータスを示します。



シリアル番号

図 10 ステータスインジケータの位置

電源インジケータ

電源インジケーターは、主電源スイッチに組み込まれています。このイン ジケーターが点灯(緑)しているときは、電源が**オン**になっています。

ステータスインジケータ

モジュールステータスインジケータ

モジュールのステータスインジケータは、次の 6 つの起こり得るモジュー ル状態の 1 つを示します。

- ステータスインジケータがオフ (電源ランプはオン)の場合は、モジュールはプレラン状態になており、分析を開始する準備が完了しています。
- 緑色のステータスインジケータは、モジュールが分析を実行中であることを示します(ランモード)。
- 黄色のインジケータは、ノットレディ状態を示します。指定状態への到 達または指定状態への完了を待機しているとき (例えば、設定値を変 更した直後)、またはセルフテスト手順の実行中は、モジュールはノッ トレディ状態になります。
- エラー状態を示す時は、ステータスインジケータは赤になります。エラー状態は、モジュールの正常な動作に影響を与える内部の問題が検出されたことを示します。通常、エラー状態には注意が必要です(リーク、内部コンポーネントの故障など)。エラー状態が発生すると、分析は中断されます。

解析中にエラーが発生すると、LC システム内に通知されるため、赤色 LED が別のモジュールの問題を示すことがあります。ユーザーインター フェイスのステータス表示を使えば、エラーの主要因 / モジュールが分 かります。

- ・ 点滅インジケータは、モジュールがレジデントモード(メインファーム ウェアの更新中など)であることを示します。
- 高速点滅インジケータは、モジュールが低レベルのエラーモードである ことを示します。このような場合は、モジュールを再起動するか、コー ルドスタートを行ってみてください(『「特別な設定」247 ページ 図』を 参照)。その後、ファームウェア更新を試します(『「モジュールファー ムウェアの交換」185 ページ 図』を参照)。問題が解決しない場合は、 メインボードの交換が必要です。

使用可能なテストとユーザーインタフェース

使用可能なテストとユーザーインタフェース

- ユーザーインターフェイスに応じて、使用できるテストとスクリーン / レポートは変わる可能性があります(「テスト機能とキャリブレーション」を参照してください)。
- 最適なツールは Agilent Lab Advisor ソフトウェアです (『「Agilent Lab Advisor ソフトウェア」82ページ 図』を参照)。
- Agilent ChemStation B. 04.02 以降には、メンテナンス / テスト機能が含まれない場合があります。
- これらの手順で使用されるスクリーンショットは Agilent Lab Advisor ソフトウェアに基づいています。

Agilent Lab Advisor ソフトウェア

Agilent Lab Advisor ソフトウェア

Agilent Lab Advisor ソフトウェアは、データシステムとは別に使用でき るスタンドアローン製品です。Agilent Lab Advisor ソフトウェアは、高 品質のクロマトグラフ結果を得るためのラボ管理に役立ち、1 台の Agilent LC、またはラボのイントラネットに設定されたすべての Agilent LC をリアルタイムでモニタリングできます。

Agilent Lab Advisor ソフトウェアは、すべての Agilent 1200 Infinity シリーズのモジュールに対する診断能力があります。これには、すべての メンテナンスルーチンに対する診断機能、キャリブレーション手順、メン テナンスルーチンが含まれます。

Agilent Lab Advisor ソフトウェアにより、ユーザーは LC 機器のステー タスをモニタリングすることもできます。EMF (Early Maintenance Feedback)機能は、予防メンテナンスの実施に役立ちます。さらに、ユー ザーは各 LC 機器のステータスレポートを作成できます。Agilent Lab Advisor ソフトウェアで提供されるテストや診断機能は、このマニュアル の説明と異なる場合があります。詳細は、Agilent Lab Advisor ソフト ウェアのヘルプファイルを参照してください。

Lab Advisor Basic は Lab Advisor ソフトウェアの基本機能バージョン で、設置、使用、メンテナンスに必要な機能は限定されています。修理、 トラブルシューティング、モニタリングなどの高度な機能は含まれていま せん。



Agilent 1290 Infinity バイナリポンプ ユーザーマニュアル

, エラー情報

7

エラーメッセージ内容 85 一般エラーメッセージ 86 Timeout 86 Shutdown 87 Remote Timeout 88 Lost CAN Partner 89 Leak Sensor Short 90 91 Leak Sensor Open Compensation Sensor Open 92 Compensation Sensor Short 92 Fan Failed 93 Leak 94 Open Cover 94 Cover Violation 95 ポンプのエラーメッセージ 96 Pressure of binary pump above upper limit 96 Pressure below lower limit 97 Target pressure not reached for binary pump degasser 98 Degasser's pressure limit violation 98 Solvent counter exceeded limit 99 Waste counter limit exceeded 100 Flow rate limit exceeded 101 Binary pump shutdown during analysis 102 Reading the pump encoder tag failed 102 Writing the pump encoder tag failed 103 Pump drive blocked or encoder failed 103 Drive current too low 104



7 エラー情報

Agilent Lab Advisor ソフトウェア

Drive Encoder failed 104 Drive current too high 105 Drive timeout 105 Overcurrent of pump drive 106 Overcurrent of solvent selection valve (SSV) 106 Deliver underrun 107 Defect connection between main board and pump drive encoder 107 Pump drive encoder defect 108 Purge valve failed 108 Reading of purge valve tag failed 109 Pump drive encoder rollover 109 Drive position limit 110 Insufficient power of drive encoder LED 110 Drive encoder error 111 Writing the purge valve tag failed 111 Current of primary pump drive too high 112 Current of secondary pump drive too high 113 Unknown purge valve type 114 Pump drive encoder error 114 Pump drive error 115 Pump drive stroke blocked 115 Pump drive stop not found 116 Pressure sensor calibration wrong or missing 116 Seal wash pump was missing when tried to turn 117 on

この章では、エラーメッセージの意味を解説し、考えられる原因に 関する情報とエラー状態から回復するための推奨方法を示します。

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。



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。

Timeout

Error ID: 0062

タイムアウト

タイムアウト値を超えました。

1 分析が正常終了した後、要求通り

にタイムアウト機能によってモ

2 シーケンスまたはマルチ注入測定

中に、タイムアウト値より長い時

間、ノットレディ状態が続いた。

ジュールをオフにしました。

考えられる原因

対策

ログブックを確認して、ノットレ 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 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。

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。



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)。

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。



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。

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

対策

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。

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。

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

対策

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。

Open Cover

Error ID: 0205

カバーが開いています

上部発泡材が取り外されました。

考えられる原因

- 対策
- 発泡材によってセンサーが有効に Agilent Technologies に連絡してくなっていません。
 ださい。
- センサーが汚れているか、故障し Agilent Technologies に連絡してくている。

Cover Violation

Error ID: 7461

カバー違反

上部発泡材が取り外されました。

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。ランプ点灯中に発泡材が取り外されると(あるいは、発泡材が取 り外された等の状態でランプの点灯を試みると)、ランプは消灯し、エ ラーメッセージが生成されます。

考えられる原因

- 1 操作中に上部発泡材が取り外され Agilent のサービス担当者に連絡し ました。 てください。
- 発泡材によってセンサーが有効に Agilent のサービス担当者に連絡しなっていません。
 なっていません。

7 エラー情報 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です。

Pressure of binary pump above upper limit

Error ID: 22014

バイナリポンプの圧力が上限値を超える

- 圧力が圧力上限値を超えました。
- パラメータ:実測圧力

考えられる原因

ている。

対策

- 1 圧力センサーの後の流路が詰まっ ・ LC システム内の詰まりをチェッ クしてください。例えば、パージ バルブ、ジェットウィーバー、劣 化したカラム、カラムフリット、 ニードル、ニードルシート、キャ ピラリなどです。
 - 溶媒内の粒子をチェックしてくだ さい。
- **2** 設定(圧力限界値、流量など)が ・ 流速を下げてください。 不適切。
 ・ 圧力リミットを上げてください。

エラー情報 7 プのエラーメッセージ

ポンプのエラーメッセージ

Pressure below lower limit

Error ID: 22015

圧力が下限値を下回った

圧力が圧力下限値を下回りました。

パラメータ:なし

考えられる原因

- **1** リーク
- 2 ボトルが空になった
- 3 間違った溶媒(粘度)の使用

4 設定が不適切

5 カラムの劣化

対策

- リークをチェックしてください。
- 溶媒ボトル容量をチェックしてくだ さい。
- 溶媒をチェックしてください。

流速と低圧力リミットをチェックし てください。

カラムを交換してください。



Target pressure not reached for binary pump degasser

Error ID: 22031

バイナリポンプのデガッサのターゲット圧力に達していない

期待される RT 内に、バイナリポンプのデガッサのターゲット圧力に達し ていません。

• パラメータ: 圧力 mbar

考えられる原因

対策

- 温度の変動により、デガッサチャ モジュールを平衡化し再起動してく ンバ内に気泡発生。
 ださい。
- **2** デガッサの故障。 Agilent のサービス担当者に連絡し てください。

Degasser's pressure limit violation

Error ID: 22032

デガッサの圧カリミット違反

圧力が限度を大きく超過

考えられる原因

- デガッサチャンバまたはデガッサ Agilent のサービス担当者に連絡し チューブのリーク。
 てください。
- 2 真空ポンプの不具合。
 Agilent のサービス担当者に連絡してください。
- デガッサチャンバが空、またはそ 未使用のデガッサチャンネルをブの接続先が大気圧。
 ロックしてください。

エラー情報 7 ポンプのエラーメッセージ

Solvent counter exceeded limit

Error ID: 22055

溶媒カウンタが制限を超えた

溶媒容量のカウンタが、ユーザー インターフェイスで設定された制限を超 えました。

パラメータ:

- ・ 溶媒選択バルブなし:
 - 0 (チャンネル A)、1 (チャンネル B)
- 溶媒選択バルブあり:

2 (チャンネル A1)、3 (チャンネル B1)、4 (チャンネル A2)、5 (チャ ンネル B2)

考えられる原因

対策

- **1** 溶媒がない。 溶媒ボトルを補充してください。
- 2 設定が不適切。 ユーザー インターフェイスの溶媒カ ウンタ設定をチェックしてください。



Waste counter limit exceeded

Error ID: 22056

廃液カウンタの制限を超えた

廃液容量のカウンタが、ユーザー インターフェイスで設定された制限を超 えました。

・ パラメータ:なし

考えられる原因

- 1 廃液コンテナがいっぱい。
- 2 廃液コンテナの設定が不適当。

対策

廃液コンテナを空にしてください。

- 廃液カウンタをリセットしてください。
- 廃液カウンタの制限を調節してく ださい。

エラー情報 7 ポンプのエラーメッセージ

Flow rate limit exceeded

Error ID: 22064

流速制限を超えた

圧力テスト等のポンプが圧力コントロールモードで動作している際に、バ イナリポンプの流量が制限を超えました。

・ パラメータ:なし

考えられる原因

対策

- リーク ポンプと流路内のリークをチェック してください。
 ボトルが空です。 溶媒ボトルを補充してください。
 シャットオフバルブが閉じていま シャットオフバルブを開いでくださ
- シャットオノハルフか闭していま シャットオノハルフを開いでくたさす。
- 4 圧力センサーでドリフトが発生 圧力センサーを交換してください。
 (所要時間が数分間の短いテストでは稀)。



Binary pump shutdown during analysis

Error ID: 22065

バイナリポンプが解析中にシャットダウン

バイナリポンプが解析中にコントロールソフトウェアまたはコントロール モジュールによりシャットダウンしました。

パラメータ:0 はオフ、1 はスタンバイ。

考えられる原因

対策

1 ポンプがシャットダウンされた。 ポンプを再起動してください。

Reading the pump encoder tag failed

Error ID: 22402

ポンプエンコーダタグの読み取りが失敗した

ポンプエンコーダタグの読み取りが失敗しました。

パラメータ:1-4 ポンプドライブについて

考えられる原因

- エンコーダとメインボードの間の Agilent のサービス担当者に連絡し 接続不良。
 なください。
- タグが見つからないか間違ってい Agilent のサービス担当者に連絡しる。または、タグとエンコーダの てください。 間の接続不良。

Writing the pump encoder tag failed

Error ID: 22405

ポンプエンコーダタグの書き込みが失敗した

ポンプエンコーダタグの書き込みが失敗しました。

パラメータ:1-4 ポンプドライブについて

考えられる原因

対策

- エンコーダとメインボードの間の Agilent のサービス担当者に連絡し 接続不良。
 てください。
- タグが間違っています。または、 Agilent のサービス担当者に連絡し タグとエンコーダの間の接続不良 てください。 です。

Pump drive blocked or encoder failed

Error ID: 22406

ポンプドライブが詰まったかまたはエンコーダが機能しなくなった

ポンプドライブが詰まったかまたはエンコーダが機能しなくなりました ・ パラメータ:なし

考えられる原因

対策

 ポンプドライブが詰まっている。 Agilent のサービス担当者に連絡し または、ドライブエンコーダの故 てください。 障。



Drive current too low

Error ID: 22407

ドライブの電流が少なすぎる

ポンプドライブの電流消費が少なすぎます。

パラメータ:1-4 ポンプドライブについて

考えられる原因

対策

- **1** ドライブモータの故障。 Agilent のサービス担当者に連絡し てください。
- ポンプドライブからメインボード Agilent のサービス担当者に連絡し への接続が間違っているか、接続 てください。 されていない。

Drive Encoder failed

Error ID: 22408

ドライブエンコーダが機能しなくなった

ドライブエンコーダが、ポンプドライブのキャリブレーション中に機能しなくなりました。

考えられる原因

1 内部のエラー。

対策

Agilent サポートに連絡してください。

エラー情報 7 ポンプのエラーメッセージ

Drive current too high

Error ID: 22409

ドライブの電流が高すぎる

ポンプドライブの電流消費が高すぎます。

パラメータ:1-4 ポンプドライブについて

考えられる原因

対策

 圧力センサの前でシステムが詰 まっています。
 アウトレットバルブ フィルタフリッ ト、パージバルブ、熱交換器などの 詰まりをチェックしてください。
 ドライブモータの故障。
 Agilent のサービス担当者に連絡し てください。

Drive timeout

Error ID: 22410

ドライブタイムアウト

ドライブが機械的に詰まると、初期化中に機能しなくなります。

パラメータ:1-4 ポンプドライブについて

考えられる原因

対策

 ポンプドライブが詰まっている。 Agilent のサービス担当者に連絡し またはドライブモータの故障。 てください。



Overcurrent of pump drive

Error ID: 22411

ポンプドライブの過電流

ポンプドライブの電流消費が高すぎます。

パラメータ:1-4 ポンプドライブについて

考えられる原因

対策

 圧力センサの前でシステムが詰 まっています。
 アウトレットバルブ フィルタフリット、パージバルブ、熱交換器などの 詰まりをチェックしてください。
 ドライブモータの故障。
 Agilent のサービス担当者に連絡してください。

Overcurrent of solvent selection valve (SSV)

Error ID: 22412

溶媒切り替えバルブ(SSV)の過電流

溶媒切り替えバルブ(SSV)の過電流。

パラメータ:なし

考えられる原因

対策

1 バルブが故障しています。 溶媒切り替えバルブの交換。

エラー情報 7 ポンプのエラーメッセージ

Deliver underrun

Error ID: 22413

送液のアンダーラン

- 内部のエラー。
- ・ パラメータ:なし

考えられる原因

1 内部のエラー。

対策

Agilent のサービス担当者に連絡し てください。

Defect connection between main board and pump drive encoder

Error ID: 22414

メインボードとポンプドライブエンコーダの接続不具合

メインボードとポンプドライブエンコーダの接続不具合

パラメータ:1-4 ポンプドライブについて

考えられる原因

- メインボードとポンプドライブエ Agilent のサービス担当者に連絡し ンコーダの接続不良。
 てください。
- **2** エンコーダの故障。 Agilent のサービス担当者に連絡してください。



Pump drive encoder defect

Error ID: 22415

ポンプドライブエンコーダの不具合

ポンプドライブエンコーダの不具合

パラメータ:1-4 ポンプドライブについて

考えられる原因

対策

1 エンコーダの故障。

Agilent のサービス担当者に連絡し てください。

Purge valve failed

Error ID: 22417

パージバルブの故障

パージバルブエンコーダの手順喪失。

パラメータ:なし

考えられる原因

- ブロックされているか、故障して をチェックしてください。 います。
- 1 パージバルブドライブが機械的に ・ パージバルブのヘッドの取り付け
 - Agilent のサービス担当者に連絡 してください。
Reading of purge valve tag failed

Error ID: 22420

パージバルブのタグの読み取り失敗

パージバルブのタグの読み取りに失敗しました。

パラメータ:なし

考えられる原因

対策

- パージバルブのタグの読み取りに ケーブルの接続をチェックしてくだ 失敗しました。
- パージバルブヘッドのタグが間 パージ バルブのヘッドを交換してく 違っているか空です。
 ださい。
- **3** パージバルブタグリーダーが故障 Agilent のサービス担当者に連絡ししています。 てください。

Pump drive encoder rollover

Error ID: 22424

ポンプドライブエンコーダのロールオーバー

無効のポンプドライブエンコーダが検出されました。

パラメータ:1-4 ポンプドライブについて

考えられる原因

対策

 ポンプドライブエンコーダの故 Agilent のサービス担当者に連絡し 障。

109



Drive position limit

Error ID: 22425

ドライブポジションリミット

内部のエラー。

パラメータ:1-4 ポンプドライブについて

考えられる原因

対策

1 内部のエラー。

Agilent のサービス担当者に連絡し てください。

Insufficient power of drive encoder LED

Error ID: 22426

ドライブエンコーダ LED の電力が不十分

ドライブエンコーダ LED の電力が不十分

パラメータ:1-4 ポンプドライブについて

考えられる原因

対策

 ポンプドライブエンコーダの故 Agilent のサービス担当者に連絡し 障。

エラー情報 7 ポンプのエラーメッセージ

Drive encoder error

Error ID: 22427- 22430

ドライブエンコーダエラー

ポンプドライブエンコーダにエラーが発生しました。

パラメータ:1-4 ポンプドライブについて

考えられる原因

対策

 ポンプドライブエンコーダの故 Agilent のサービス担当者に連絡し 障。
 Cください。

Writing the purge valve tag failed

Error ID: 22431

パージバルブのタグの書き込み失敗

パージバルブのタグの書き込みが失敗しました。

・ パラメータ:なし

考えられる原因

対策

- パージバルブヘッドのタグが間 パージ バルブのヘッドを交換してく 違っています。
 ださい。
- パージバルブタグリーダーが故障 Agilent のサービス担当者に連絡ししています。



Current of primary pump drive too high

Error ID: 22433

プライマリポンプドライブの電流が高過ぎます

プライマリポンプドライブの電流が高すぎます。

考えられる原因

対策

- 1 熱交換器など、プライマリポンプ・流路の詰まりを確認してくださ ヘッドと圧力センサの間の流路が 詰まっています。
- しています。
- い。
- Agilent のサービス担当者に連絡 してください。
- 2 プライマリポンプドライブが故障 Agilent のサービス担当者に連絡し てください。

Current of secondary pump drive too high

Error ID: 22434

セカンダリポンプドライブの電流が高過ぎます

セカンダリポンプドライブの電流が高すぎます。

パラメータ:2 または3 ポンプドライブについて。

考えられる原因

対策

- 1 熱交換器など、セカンダリポンプ・ 流路の詰まりを確認してくださ ヘッドと圧力センサの間の流路が 詰まっています。
- しています。
- い。
- Agilent のサービス担当者に連絡 してください。
- 2 セカンダリポンプドライブが故障 Agilent のサービス担当者に連絡し てください。



Unknown purge valve type

Error ID: 22435

不明なパージバルブ タイプ

- パージバルブのタイプ情報が無効です。
- パラメータ:なし

考えられる原因

対策

るか、交換してください。

- **1** 間違ったバルブヘッドが取り付け パージバルブのヘッドをチェックす られています。
- 2 バルブヘッドの RFID タグの内容 パージバルブのヘッドをチェックす るか、交換してください。 が無効です。

Pump drive encoder error

Error ID: 22437

ポンプドライブエンコーダのエラー

ポンプドライブエンコーダがシグナルを作成しませんでした。

パラメータ:1-4 ポンプドライブについて

考えられる原因

対策

1 ポンプドライブエンコーダの故 Agilent のサービス担当者に連絡し てください。 障。

エラー情報 7 ポンプのエラーメッセージ

Pump drive error

Error ID: 22438, 22439

ポンプドライブのエラー

ポンプドライブが、キャリブレーション中に機能しなくなりました。

パラメータ:1-4 ポンプドライブについて

考えられる原因

対策

ポンプドライブモータが故障して Agilent のサービス担当者に連絡しいるか、機械的にブロックされて てください。います。

Pump drive stroke blocked

Error ID: 22441

ポンプドライブストロークがブロックされた

ポンプドライブの動作がブロックされました。

パラメータ:1-4 ポンプドライブについて

考えられる原因

対策

 ポンプドライブモータが機械的に Agilent のサービス担当者に連絡し ブロックされている。
 てください。



Pump drive stop not found

Error ID: 22442

ポンプドライブストップが見つからない

ポンプドライブストップが見つかりません。

パラメータ:1-4 ポンプドライブについて

考えられる原因

対策

1 ポンプドライブスピンドルの故 Agilent のサービス担当者に連絡し てください。 障。

Pressure sensor calibration wrong or missing

Error ID: 22443

圧力センサのキャリブレーションが誤っているか、センサがありません

圧力センサのキャリブレーションが誤っているか、センサがありません。 パラメータ:なし

考えられる原因

対策

- 1 圧力センサのキャリブレーション ・ 圧力センサーを交換してくださ が間違っているか、行われていま せん。
 - V.
 - Agilent のサービス担当者に連絡 してください。

Seal wash pump was missing when tried to turn on

Error ID: 22499

オンにしようとしてもウォッシュ用シールポンプがない

ウォッシュ用シールポンプが検出されません(以前は設定されたり、また は検出されましたが)

考えられる原因

対策

- ウォッシュ用シールポンプへのケーブルの接続をチェックしてくだケーブル接続の不具合。
- 2 ウォッシュ用シールポンプモー Agilent のサービス担当者に連絡し タの不具合。
- **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。

エラー情報 ポンプのエラーメッセージ

Agilent 1290 Infinity バイナリポンプ ユーザーマニュアル



8 テスト機能とキャリブレーション

はじめに 120 システム圧力テスト 121 システム圧力テストの評価 123 ポンプヘッドのリークテスト 124 ポンプヘッドのリークテストの評価 126

この章では、モジュールのテストについて説明します。



8 テスト機能とキャリブレーション はじめに

はじめに

すべてのテストの説明は、Agilent Lab Advisor ソフトウェア B.01.04 に 基づくものです。他のユーザーインタフェースについては、まったくテス トされていないか、あるいは部分的にしかテストされていません。

インタフェース コメント 使用可能な機能 アジレント機器ユー メンテナンステストを システム圧力テスト ティリティ 提供 Agilent Lab Advisor すべてのテストを提供 システム圧力テスト ポンプヘッドのリー クテスト Agilent ChemStation 提供されるテストなし 圧力 クロマトグラフ信号へ の圧力の追加が可能 温度メインボード Agilent Instant Pilot 一部のテストを提供 システム圧力テスト 値の監視
 E力
 • 圧力リップル 流量(動作圧力の) 場合)

表 8 インタフェースと使用可能なテスト機能

インタフェースの使用の詳細については、そのインタフェースのマニュア ルを参照してください。

テスト機能とキャリブレーション 8

システム圧力テスト

システム圧力テスト

このテストでは、ポンプのアウトレットバルブとブランクナットの間でシ ステムの漏れ率を測定します。システム内のフローセルよりも前のさまざ まな位置にブランクナットを配置することで、個々のモジュールとコン ポーネントの漏れ率を測定することができます。このテストでは、テスト を実施する圧力を設定できます。高圧部品の漏れ率は必ずしも線形関数に ならないため、システムの通常の動作圧力に対応する圧力でテストを実施 することをお勧めします。

- **日時**: リークが疑われる場合。メンテナンスタスクが正しく実行されたことを確認 するため。
- **必要な部品: 番号 部品番号 説明** 1 01080-83202 ブランクナット
- **必要な準備:** 両方のチャンネルに溶媒が含まれている必要があります。

- 8 テスト機能とキャリブレーション システム圧力テスト
 - 推奨するユーザーインタフェースを使用して System Pressure Test 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へ ルプを参照)。

Test Name Module Approx Time		ime Time	System Pressure Test G4220A:LP00000005 5 min	Description No description		
Sta	Status		Running			
Test	Proc	edure —		Result		
	1	Proporo	nump pressure test	Name	Value	
	1.	r iepaie	pump pressure rest	Maximum system pressure	1200 bar	
	2.	Enter the	e test pressure	Channel A2	Done	
\checkmark	3.	Configur	re purge process and parameter	Remaining purge time	0 min 0 sec	
1	4.	Insert bl	ank nut			
u))	5.	System	checking leak rate of system			
	6.	Evaluate	e results			
	7.	Restore	system configuration			

図 11 システム圧力テスト - 結果



テスト機能とキャリブレーション 8

システム圧力テスト

システム圧力テストの評価

Test Failed

テスト失敗

対策 考えられる原因 **1** ブランクナットの損傷 (締めすぎ 考えられる他の失敗の原因を調べる 前に、使用しているブランクナット による変形) の状態が正常で、適切に締められて いることを確認してください。 ポンプヘッドのリークテストを実施 2 ポンプのリーク します。 3 フィッティングの緩みまたはリー フィッティングを締めるか、キャピ ラリを交換します。 ク 4 オートサンプラのリーク オートサンプラのリークテストを実 施します。 5 カラムコンパートメントのバルブ TCC バルブのロータシールを交換し ます。 のリーク

注記 テストのエラーと結果の失敗との違いに注意してください。エラーは、テスト 実行中の異常終了により発生しますが、結果の失敗は、テスト結果が指定され た限界値内になかったことを示します。

8 テスト機能とキャリブレーション ポンプヘッドのリークテスト

ポンプヘッドのリークテスト

このテストでは、各ポンプヘッドを個別にブロックし、指定レベルまで加 圧することにより、個々のポンプヘッドのリークを確認します。フローが セカンダリピストンによって提供され、アウトレットボールバルブとパー ジバルブの間の部品の耐圧性がテストされます。このテストでは、テスト を実施する圧力を設定できます。高圧部品の漏れ率は必ずしも線形関数に ならないため、システムの通常の動作圧力に対応する圧力でテストを実施 することをお勧めします。

日時: 過度の圧力リップルがある場合や、ポンプ性能の問題が疑われる場合。

必要な準備: 両方のチャンネルに溶媒が含まれている必要があります。

テスト機能とキャリブレーション 8

ポンプヘッドのリークテスト

 推奨するユーザーインタフェースを使用して Pump Head Leak Test 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)。

	st Na	me	Pump Head Leak Test	Description No description			
Module			G4220A:LP00000005				
Ap	Approx. Time Status		5 min Running				
Sta							
est	: Proc	edure —		Result Name	Value		
est	: Proc	edure —					
	1.	Checkir	ng pre-requisites (pump on etc.)	Name Maximum system pressure	1200 bar		
	2.	Enter th	e test pressure	Channel A2	Done		
	З.	Configu	re purge process and parameter	Channel B2	Done		
1	4.	System	checking leak tightness in channel A	Remaining purge time	0 min 0 sec		
•		-	obooking look tightnoon in obonnol P				
¢	5.	System	checking leak ughtness in channel b				
\$	5. 6.	System Evaluat	e results				



8 テスト機能とキャリブレーション ポンプヘッドのリークテスト

ポンプヘッドのリークテストの評価

Test Failed

テスト失敗

考えられる原因

対策

- フィッティングの緩みまたはリーク フィッティングを締めるか、キャピ ラリを交換します。
- 2 高圧フィルタアセンブリのリーク ・ フィルタのハウジングとフィッ
 - ティングを締めます。
 - 注 テストに成功する可能性を 上げるために、加圧していないシ ステムのフィッティングだけを締 めてください。
 - テストを再び実施します。
- ポンプシールまたはピストンの損傷
 ポンプヘッドまたはポンプシール を交換します。
 - 注 ポンプシールが間違って取り付けられていると、ポンプシールの寿命が大幅に短縮します。
- 4 アウトレットボールバルブのリーク ポンプヘッドまたはアウトレット ボールバルブを交換します。
- 5 パージバルブのリーク パージバルブのロータシールを交換 します。

テストの**エラー**と結果の失敗との違いに注意してください。エラーは、テスト 実行中の異常終了により発生しますが、結果の失敗は、テスト結果が指定され た限界値内になかったことを示します。

注記



Agilent 1290 Infinity バイナリポンプ ユーザーマニュアル

-メンテナンス

9

メンテナンスの概要 128 警告と注意 130 メンテナンスの概要 132 モジュールのクリーニング 133 フィッティングとキャピラリの設置 134 シャットオフバルブパネルの交換 135 圧力センサーの交換 137 インレットバルブの交換 139 アウトレットバルブの交換 141 溶媒切り替えバルブ(SSV)の交換 144 構成の変更または Jet Weaver の交換 146 シールウォッシュポンプの交換 149 スタックインレットバルブの取り外し 150 ポンプヘッドの交換 153 ポンプヘッドの分解 159 プライマリポンプヘッドの分解 161 セカンダリポンプヘッドの分解 164 ポンプヘッドの組み立て 167 パージバルブヘッドの交換 177 高圧フィルタアセンブリの部品の交換 180 バルブレールキットの設置 182 主電源ヒューズの交換 183 モジュールファームウェアの交換 185 ポンプモジュールの移動の準備 186

この章では、Agilent 1290 Infinity バイナリポンプのメンテナン スについて説明します。



9 メンテナンス メンテナンスの概要

メンテナンスの概要

『128 ページ 図 15』メインユーザーがアクセスできる Agilent 1290 Infinity バイナリポンプのアセンブリを示します。これらの部品は前面から簡単に修理できます。システムスタックからポンプを取り外す必要はありません。



パージバルブ

図 15 メンテナンス部品

『129ページ図16』に、これらのメインアセンブリ間の配管を示します。

メンテナンス 9 メンテナンスの概要



9 メンテナンス 警告と注意

警告と注意

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。

警告 感電

モジュールの修理作業によって人身障害が起こる恐れがあります(カバーを開けたままにして感電するなど)。

- → 本装置のカバーは取り外さないでください。
- → モジュール内部の修理は、有資格者だけに許可されています。

警告

人身障害と製品の損害

アジレントは、全部または一部において、製品を不正に利用したり、 製品を許可なく改変、調整、修正した場合、アジレント製品ユー ザーガイドに従わなかった場合、または適用される法律、法令に違 反して製品を使用した場合に生じるいかなる損害にも責任を負いま せん。

→ アジレント製品は、アジレント製品ユーザーガイドに記載された 方法で使用してください。

メンテナンス 9 警告と注意

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。

メンテナンスの概要

メンテナンスの概要

以降のページでは、メインカバーを開けずに実施できるモジュールのメン テナンス (簡単な修理) について説明します。

メンテナンス 9 モジュールのクリーニング

モジュールのクリーニング

モジュールケースを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。

警告
 モジュールの電子コンパートメントに液体が入ると、感電やモジュールの損傷を引き起こす恐れがあります。
 → 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。

警告

フィッティングとキャピラリの設置

フィッティングとキャピラリの設置

| 溶媒は高圧下で噴出できます。

- → 流路を開くときは、適切な安全手順を守ってください(例えば、 ゴーグル、安全手袋、防護衣着用)。
- - → 圧力下では決して配管を締めないでください。
 - 注記 フィッティングの寿命は、それが締められている固さ次第です。固く締められ ていると寿命は縮まります。
 - フィッティングを締めすぎた場合、交換してください。
 - 1 フィッティングとキャピラリを取り付けます。
 - **2** フィッティングとキャピラリを締めます。

メンテナンス 9 シャットオフバルブパネルの交換

シャットオフバルブパネルの交換

- **日時**: シャットオフバルブが破損した場合、または修理手順のためにパネルを取り 外す必要がある場合。
- 必要な部品: 番号 部品番号 説明
 1 5067-4124 シャットオフバルブ
 1 G4220-40004 シャットオフバルブパネル
 2 G4220-60035 チューブキット140 mm SSV からシャットオフバルブ またはデガッサユニットまで(チューブ2本)

必要な準備: リークを防ぐために、チューブを溶媒ボトルから取り外します。



シャットオフバルブパネルの交換



メンテナンス 9 圧力センサーの交換

圧力センサーの交換

日時: 圧力シグナルがない、または無効

必要なツール: 部品番号 説明 8710-2412 六角レンチ 2.5 mm、15 cm 長、ストレートハンドル付き 8710-0510 1/4 インチ × 5/16 インチのレンチ スクリュードライバー

必要な部品: 番号 部品番号 説明 1 G4220-60001 圧力センサー 120 MPa (1200 bar)

必要な準備: ポンプ流量をオフにし、ポンプの電源を切ります。

注記 ここでは、圧力センサーの交換手順を説明します。 センサーへのに接続するケーブルの交換も必要な場合は、Agilent のサービス 担当者に連絡してください。



圧力センサーの交換



インレットバルブの交換

インレットバルブの交換

日時: インレットバルブが故障している場合。

必要なツール: 部品番号 説明 レンチ、14 mm G4220-20012 トルクレンチ 2 - 25 Nm

必要な部品: 部品番号 説明 G4220-60022 インレットバルブ (プライマリポンプヘッド)

注意 インレットバルブの潜在的な損傷
 きつく締めすぎるとバルブを損傷し、トルクがゆるいと漏れます。
 → インレットバルブの修理には、トルクレンチ(10 Nm)を使用します。



インレットバルブの交換



アウトレットバルブの交換

アウトレットバルブの交換

日時: アウトレットバルブが故障している場合。

必要なツール: 部品番号 説明 レンチ、14 mm G4220-20012 トルクレンチ 2 - 25 Nm G4220-20014 2.5 mm 六角ビット

必要な部品: 部品番号 説明

G4220-60028 アウトレットバルブ 1290 Infinity クォータナリポンプ (プライマリポンプヘッド)

G4220-20020 1290 Infinity アウトレットバルブ用の内部ゴールドシール



アウトレットバルブの交換



アウトレットバルブの交換



溶媒切り替えバルブ(SSV)の交換

溶媒切り替えバルブ(SSV)の交換

日時:

溶媒切り替えバルブに問題が発生した場合。

必要な部品: 番号 部品番号 説明

1 G4280-60029 溶媒切り替えバルブ

4 G4220-60035 チューブキット 140 mm SSV からシャットオフバルブ またはデガッサユニットまで(チューブ 2本)


溶媒切り替えバルブ(SSV)の交換



構成の変更または Jet Weaver の交換

構成の変更または Jet Weaver の交換

- **日時**: ポンプ設定をミキシング性能または低ディレイ ボリューム / 高速グラジエントに最適化するには、性能の最適化の章を参照してください。
- **必要なツール: 部品番号 説明** 8710-0510 1/4 インチスパナ 3 mm 六角レンチ
- 必要な部品: 番号 部品番号 説明
 1 G4220-60006 Jet Weaver 35 µL/100 µL
 1 G4220-60012 Jet Weaver 380 µL (オプション)
 1 G4220-87000 キャピラリ、SST バルブから Jet Weaver まで 300 mm x0.17 mm 内径

構成の変更または Jet Weaver の交換



構成の変更または Jet Weaver の交換



メンテナンス 9 シールウォッシュポンプの交換

シールウォッシュポンプの交換

日時: ウォッシュ用シールポンプが摩耗している場合。

必要な部品: 部品番号 説明 5042-8507 ウォッシュ用シール用ペリスタルポンプ

必要な準備: シャットオフバルブパネル(『「シャットオフバルブパネルの交換」135 ページ図』)と配管をシールウォッシュポンプから取り外します。



注記

スタックインレットバルブの取り外し

スタックインレットバルブの取り外し

日時: 長期間電源を切っておいた後にインレットバルブが詰まった場合、またはポンプが加圧できない場合。

システムの電源を長期間切断する前に、少なくとも10%の IPA でフラッシュ し、インレットバルブの詰まりを防止します。



スタックインレットバルブの取り外し



スタックインレットバルブの取り外し



メンテナンス 9 ポンプヘッドの交換

ポンプヘッドの交換

Agilent 1290 Infinity バイナリポンプには、2 つの溶媒チャンネル A と B に対して 2 つのポンプアセンブリがあり、どちらも 2 つのポンプドラ イブとポンプヘッドで構成されています。プライマリーポンプヘッドから 各ポンプに入った溶媒は、セカンダリポンプヘッドに移送され、ポンプバ ルブに接続されたセカンダリポンプヘッドのアウトレットを出ます。



図 17 ポンプヘッドアセンブリ(左)

注意

注意

ポンプヘッドの交換

日時: ポンプの性能に問題が発生した場合。

- 必要なツール:部品番号説明G4220-20012トルクレンチ 2 25 NmG4220-200134 mm 六角ビットG4220-20015アダプタ ¼ インチ、正方形 / 六角形
- 必要な部品: 番号 部品番号 説明 ポンプヘッドアセンブリチャンネル A (左) (シール G4220-60200 1 ウォッシュオプション付き) ポンプヘッドアセンブリチャンネル B (右) (シール 1 G4220-60210 ウォッシュオプション付き) ポンプヘッドアセンブリチャンネル A(左)(シール 1 G4220-60400ウォッシュオプションなし) ポンプヘッドアセンブリチャンネル B (右) (シール 1 G4220-60410 ウォッシュオプションなし)

必要な準備: ・ 主電源スイッチでポンプの電源を切ります。

• フロント カバーを取り外します。

寿命の制限

ポンプヘッドアセンブリは、標準のツールでは組み立て直しができない交換部品です。ポンプヘッドを分解すると、寿命が大きく制限されます。

→ ポンプヘッドアセンブリは分解しないでください。

接続の破損

1 つのポンプヘッドアセンブリ(溶媒チャンネル)の両方のポンプヘッド 間の配管を分解すると、接続を破損し、リークを引き起こすことがありま す。

→ ポンプヘッド間の配管は外さないでください。

メンテナンス 9 ポンプヘッドの交換



→ ポンプヘッドのネジに強い力を与えないでください。

→この目的にはトルク六角レンチを使用してください。

ポンプピストンの損傷 注意 メンテナンスポジション以外のポジションでポンプヘッドを取り外すと、 ポンプピストンが破損することがあります。 → ポンプの電源を切る前に、メンテナンスポジションに移動してくださ い。

この手順では、左のポンプヘッドアセンブリの交換を説明します。右のポンプ 注記 ヘッドアセンブリも同様に交換できます。

1 つのポンプヘッドアセンブリは 2 つのポンプヘッドから構成され、両方の ポンプヘッドは同時に取り外されます。



ポンプヘッドの交換



ポンプヘッドの交換



9 メンテナンス ポンプヘッドの交換

7 配管を取り付けます。	8 モジュールを完全に組み立て直したら、ポ
• デガッサユニットの出口をプライマリポ	ンプをパージする必要があります。10 min
ンプヘッドの入口に接続し、セカンダリ	間、10 ml/min の流量で溶媒を流します。
ポンプヘッドの出口をパージバルブの入	このときに、アプリケーションに適した混
口に接続します。	合比 50:50 の溶媒をチャンネル A および B
• チャンネル A(左側のポンプヘッドアセ	に使用します(50 % アセトニトリル、50 %
ンブリ)をポート 4 に、チャンネル B(水など)。
右側のポンプヘッドアセンブリ)をポー	
ト 1 に接続します。	

メンテナンス 9 ポンプヘッドの分解

ポンプヘッドの分解

日時: ポンプヘッド内部の部品を交換する必要がある場合。

- 必要なツール:部品番号説明G4220-20012トルクレンチ 2 25 NmG4220-200134 mm 六角ビットG4220-200142.5 mm 六角ビットG4220-20015アダプタ ¼ インチ、正方形 / 六角形8710-05101/4 インチ × 5/16 インチのレンチ
- **必要な準備:**『「ポンプヘッドの交換」153ページ図』の1~4の手順に従って、ポンプ ヘッドアセンブリを取り外します。
- 注記 この手順では、チャンネル A のポンプヘッドの交換について説明します。 チャンネル B も同様に交換できます。プライマリポンプヘッドには熱交換器 はありません。シールウォッシュ部品は、いずれのポンプヘッドの場合もオプ ションです。

ポンプヘッドの分解



メンテナンス 9 プライマリポンプヘッドの分解

プライマリポンプヘッドの分解

注意 ポンプピストンの損傷

- ポンプピストンは、ZrO₂ ベースのセラミックでできており、とても固く耐性のある材質ですが、横からのせん断力に敏感です。
- → 背面からポンプピストンを取り外さないでください。
- → ポンプ シールを取り外すためにピストンを使用しないでください。



プライマリポンプヘッドの分解



メンテナンス 9 プライマリポンプヘッドの分解

 ポンプピストンのシールを交換する際は、 ピストンに傷や溝、へこみがないか確認します。 	6 インサートツールのプラスチック側を使って、ポンプチャンバハウジングからポンプシールを取り外します。
注記 損傷のあるピストンは微細なリークを引き起こ して、シールの寿命を縮めます。	インサートツール
	ポンプシール
	注記 ポンプピストンでシールを外さないでくださ い!
7 ウォッシュ用シールを新しいものに交換す る場合は、インサートツールのスチール側 を使用して取り外してください。	
ウォッシュ用シール	



注意

セカンダリポンプヘッドの分解

📕 ポンプピストンの損傷

- ポンプピストンは、ZrO₂ ベースのセラミックでできており、とても固く耐性のある材質ですが、横からのせん断力に敏感です。
- → 背面からポンプピストンを取り外さないでください。
- → ポンプ シールを取り外すためにピストンを使用しないでください。



セカンダリポンプヘッドの分解



9 メンテナンス セカンダリポンプヘッドの分解

5 ポンプピストンのシールを交換する際は、	6 インサートツールのプラスチック側を使っ
ピストンに傷や溝、へこみがないか確認し	て、ポンプチャンバハウジングからポンプ
ます。	シールを取り外します。
注記 損傷のあるピストンは微細なリークを引き起こ して、シールの寿命を縮めます。	インサートツール
	ポンプシール
	注記 ポンプピストンでシールを外さないでくださ い!
7 ウォッシュ用シールを新しいものに交換す	
る場合は、インサートツールのスチール側 を使用して取り外してください	
ウォッシュ用シール	

ポンプヘッドの組み立て

ポンプヘッドの組み立て

日時: ポンプヘッドを取り付ける前。

- 必要なツール: 部品番号 説明 ポンプヘッドアライメントツール
 G4220-20012 トルクレンチ 2 - 25 Nm
 G4220-20013 4 mm 六角ビット
 G4220-20014 2.5 mm 六角ビット
 G4220-20015 アダプタ ¼ インチ、正方形 / 六角形
 01018-23702 インサートツール
- 必要な部品:部品番号説明0905-1420PE シール (2 個入)0905-1718ウォッシュ用シール PE5062-2484ガスケット、ウォッシュ用シール用 (6 個入)

詳細については、「部品」を参照してください。

注意 ポンプヘッド寿命の制限

バックアップシールを間違って挿入すると、ポンプヘッドの寿命が制限されることがあります。

→ バックアップシールの正しい向きに注意してください。

注意 ポンプピストンの損傷

ポンプピストンは、横からのせん断力に非常に敏感です。

→ ポンプヘッドアライメント ツールのアライメントピストンを使用して、 以下に説明する手順でアライメントを行います。

ポンプヘッドの組み立て

注意サポートリング上のピンの誤った向き

- サポートリング上のピンの正しい向きに注意せずにポンプヘッドを組み立 てると、ピストンやポンプヘッドにリークや損傷が生じることがあります。
- → サポートリング上のピンに注意してください。ポンプヘッドの部品を正しい向きに組み立てるのに役立ちます。

注意 ポンプヘッドアセンブリを取り付ける時、ポンプドライブをメンテナンス ポジションに移動させる必要があります。デフォルトポジションでポンプ ドライブを使用すると、ポンプヘッドアセンブリが損傷します。 → ポンプドライブをメンテナンスポジションにします。

注記 ここでは、ポンプヘッド アライメントツールを使用して、セカンダリポンプ ヘッドの組み立て方の手順を説明します。プライマリポンプヘッドの組み立て は適宜行うことができます。セカンダリポンプヘッドは熱交換器キャピラリを もち、これはアライメントツールにフィットしなくてはなりません。一方、プ ライマリポンプヘッドには熱交換器は含まれません。

ポンプヘッドの組み立て



ポンプヘッドの組み立て



ポンプヘッドの組み立て



ポンプヘッドの組み立て



ポンプヘッドの組み立て



ポンプヘッドの組み立て



ポンプヘッドの組み立て



ポンプヘッドの組み立て



メンテナンス 9 パージバルブヘッドの交換

パージバルブヘッドの交換

日時:

パージバルブに問題が発生した場合。

必要な部品:	番号	部品番号	説明
	1	5067-4119	パージバルブヘッド
	1	5067-4655	ポンプヘッドアセンブリチャンネル A および B への キャピラリ(2 本)
	1	G4220-87000	キャピラリ、SST バルブから Jet Weaver まで 300 mm x0.17 mm 内径
	1	5067-4656	圧力センサーへのキャピラリ(2 本)

必要な準備: パージバルブへのすべてのキャピラリ接続を取り外します。

注意

バルブヘッドを損傷したりバルブが故障したりする恐れ ポンプがオンになるとバルブタグがアクセス(読み取り / 書き込み)され、 バルブを正しく位置決めするのに使用されます。

ポンプがオンの間にバルブヘッドが交換されると、無効な情報がバルブ ヘッドに書き込まれて使用できなくなったり、位置が正しくないためにバ ルブ内の配管に不具合が生じ、部品を損傷する恐れが生じたりします。

→ ポンプの電源を切ってからパージバルブの作業を開始してください。

注意 偏った測定結果 バルブドライブには高感度の光学部品が含まれます。これらの部品が汚れ ているとバルブポートの正確な選択を妨げる可能性があり、そのため偏っ た測定結果になります。 → 光学部品は、埃やその他の汚染物質から保護してください。

パージバルブヘッドの交換



パージバルブヘッドの交換



179

高圧フィルタアセンブリの部品の交換

高圧フィルタアセンブリの部品の交換

- 日時: 高圧フィルタアセンブリの詰まりとリークを解決するため。アウトレットバルブのフィルタフリットは、システムの使用状況に応じて定期的に交換する必要があります。その他の部品は Agilent 点検 (PM) サービスの対象となります。
- **必要なツール: 部品番号 説明** G4220-20012 トルクレンチ 2 - 25 Nm G4220-20015 アダプタ ½ インチ、正方形 / 六角形 14 mm 六角ビット
- 必要な部品: 番号 部品番号 説明
 1 01018-22707 PTFE フリット(5 個)
 1 5067-4728 シールキャップ

注意 接続のリークまたは破損 プライマリポンプヘッドの出口を開くと、ポンプヘッド間の接続でリーク が発生したり、破損したりすることがあります。

→ プライマリポンプヘッドの出口は開かないでください。

注記	この手順はチャンネル A (左のポンプヘッドアセンブリ)の交換の説明で、 チャンネル B にも適用できます。どちらのケースも、メンテナンスが行われ るのはフィルタフリットを受け入れるセカンダリポンプヘッドアウトレットに だけです。

注記 PTFE フリットを交換するとき、漏れ防止にシールキャップも交換することを 検討してください。
高圧フィルタアセンブリの部品の交換



バルブレールキットの設置

バルブレールキットの設置

- **日時**: このレールは外部バルブの取り付けに必要です。
- **必要なツール: 説明** Pozidrive スクリュードライバー #1
- 必要な部品:番号部品番号説明15067-4634バルブレールキット
 - 注記 レールは、ポンプの左または右側に取り付けることができます。この手順では 左側の取り付けについて説明します。この手順は右側にも同様に適用されま す。
 - バルブレールは4本のネジでポンプカバーに固定されます。モジュール カバーの下側のネジ位置には印が付いています。最初にこれらのネジを 締め、次に上側のネジを締めます。



メンテナンス 9 主電源ヒューズの交換

主電源ヒューズの交換

- 日時:
 主電源ボタンを押しても主電源の LED が点灯しない場合(『「ステータスインジケータ」79 ページ図』を参照)。
- **必要なツール: 説明** マイナスドライバー
- **必要な部品: 番号 部品番号 説明** 1 2110-1004 ヒューズ 10 A t
- **必要な準備:** 機器の電源を切り、主電源ケーブルを外します。
- 警告
 火災の危険
 間違ったヒューズを使用すると、火災の危険があります。
 ・ 火災防止のため、必ずタイプと定格が同じ電源ヒューズとだけ交換してください。
 ・ この機器に指定されたヒューズだけを使用してください。他の ヒューズや材料の使用は禁止されています。

注記 装置の内側にはさらにヒューズがあります。 主電源ヒューズを交換しても問題が解決しない場合は、アジレントのサービス 担当者にご連絡ください。

主電源ヒューズの交換



モジュールファームウェアの交換

モジュールファームウェアの交換

日時: 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 LAN/RS-232 ファームウェア更新ツール または Agilent ラボアドバイザソフトウェア インスタントパイロット G4208A または (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、以下の操作を行います。

- 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。

Agilent 1290 Infinity バイナリポンプ ユーザーマニュアル

185

ポンプモジュールの移動の準備

ポンプモジュールの移動の準備

日時: モジュールを移動または輸送する必要がある場合。

必要な部品:	番号	部品番号	説明
	1	9301-0411	シリンジ (プラスチック製)
	1	9301-1337	シリンジアダプタ
	1	G4220-44000	保護用緩衝材

必要な準備: 両方の溶媒チャンネルをイソプロパノールでフラッシュします。

注意機械的な損傷

- → モジュールを輸送する場合は、保護用緩衝材を挿入し、機械的な損傷からモジュールを保護します。
- →保護用緩衝材をモジュールに挿入するときは、チューブやキャピラリ接続が破損しないように注意してください。

ポンプモジュールの移動の準備

1 溶媒インレットを溶媒ボトルから外します。両方の溶媒チャンネルのプ ライマリポンプヘッドの入口から溶媒チューブを外します。溶媒ボト ル、シャットオフバルブパネル、溶媒切り替えバルブ、デガッサユニッ ト、ポンプインレットの間の溶媒チューブから、シリンジを使用して液 体を取り除きます。必要に応じて溶媒切り替えバルブを交換します。



インレットプライマリポンプヘッド

ポンプモジュールの移動の準備



2 その他のモジュールおよび溶媒キャビネットへのチューブとキャピラリ 接続を取り外します。チューブプラグを取り外します。

アウトレットアクティブシールウォッシュ(オプション)

ポンプモジュールの移動の準備



3 シャットオフバルブパネルを下向きに引いて取り外します。

4 内部のチューブとキャピラリ接続はそのままにしておくこともできます。



圧力センサー

ポンプモジュールの移動の準備

- 5 他のモジュールへのケーブル接続を取り外します。モジュールをスタッ クから取り外します。
- 6 保護用緩衝材を機器の正面部分に慎重に挿入します。チューブやキャピ ラリ接続に損傷を与えないでください。



7 前面カバーを閉めてください。



8 移動や輸送の際は、モジュールとアクセサリキットを元の輸送用の箱に 入れます。

Agilent 1290 Infinity バイナリポンプ ユーザーマニュアル

10 メンテナンス用部品と器材

メンテナンス部品の概要192キャピラリ193溶媒キャビネットキット194シールウォッシュオプション195ポンプヘッドアセンブリの部品196プライマリポンプヘッドの部品198セカンダリポンプヘッドの部品202パージバルブ206カバー部品207リーク部品208ヒューズ208アクセサリキット209その他210

この章では、メンテナンス用部品について説明します。



Agilent Technologies

10 メンテナンス用部品と器材 メンテナンス部品の概要

メンテナンス部品の概要

部品番号	説明
G4280-60029	溶媒切り替えバルブ
G4220-60006	Jet Weaver 35 $\muL/100\muL$
G4220-60012	Jet Weaver 380 µL (オプション)
G4220-60001	圧力センサー 120 MPa (1200 bar)
G4220-60200	ポンプヘッドアセンブリチャンネル A(左)(シール ウォッシュオプション付き)
G4220-60210	ポンプヘッドアセンブリチャンネル B(右)(シール ウォッシュオプション付き)
G4220-60400	ポンプヘッドアセンブリチャンネル A(左)(シール ウォッシュオプションなし)
G4220-60410	ポンプヘッドアセンブリチャンネル B(右)(シール ウォッシュオプションなし)

メンテナンス用部品と器材 10 キャピラリ

キャピラリ

部品番号	説明
5067-4655	ポンプヘッドアセンブリチャンネル A および B への キャピラリ(2 本)
G4220-87000	キャピラリ、SST バルブから Jet Weaver まで 300 mm x0.17 mm 内径
5067-4656	圧力センサーへのキャピラリ(2 本)

溶媒キャビネットキット

溶媒キャビネットキット

部品番号	説明
5067-4644	溶媒キャビネットキット 1290 Infinity ポンプ 以下の部品が含まれます。
5065-9981	溶媒キャビネット 1200 Infinity、すべてのプラス チック製部品を含む)
9301-1420 (3x)	溶媒ボトル、透明
9301-1450	溶媒ボトル、褐色
5067-4124 (4x)	シャットオフバルブ
G4220-60007 (4x)	ボトルヘッドアセンブリ
G4220-60035 (4x)	チューブキット140 mm SSV からシャットオフバルブま たはデガッサユニットまで(チューブ 2 本)
G4220-40004	シャットオフバルブパネル
5042-9967	チューブ止め具 (止め具 5 本入)

シールウォッシュオプション

シールウォッシュオプション



図 18 シールウォッシュポンプ

部品番号	説明
5042-8507	シールウォッシュポンプカートリッジ
5065-9978	チューブ、内径1 mm 、外径3 mm 、シリコン、5 m、 ウォッシュ用シールオプション用

ポンプヘッドアセンブリの部品

ポンプヘッドアセンブリの部品



図 19 ポンプヘッドアセンブリの部品

メンテナンス用部品と器材 10 ポンプヘッドアセンブリの部品

品目	部品番号	説明
1	G4280-60026	高圧フィルタアセンブリ (セカンダリポンプヘッド)
2	G4220-60028	アウトレットバルブ 1290 Infinity クォータナ リポンプ (プライマリポンプヘッド)
3	G4220-60022	インレットバルブ (プライマリポンプヘッド)

10 メンテナンス用部品と器材 プライマリポンプヘッドの部品

プライマリポンプヘッドの部品

プライマリポンプヘッド、シールウォッシュ付き (バイナリポンプ)



図 20 プライマリポンプヘッド (バイナリポンプ)、 シールウォッシュ付き

メンテナンス用部品と器材 10 プライマリポンプヘッドの部品

品目	部品番号	説明
1	0515-1218	ネジ、M5、長さ 40 mm
2		ポンプチャンバハウジング (ポンプヘッドをご注文ください)
3	G4220-60022	インレットバルブ (プライマリポンプヘッド)
4	G4220-60028	アウトレットバルブ 1290 Infinity クォータナリ ポンプ (プライマリポンプヘッド)
5	G4220-20020	1290 Infinity アウトレットバルブ用の内部ゴール ドシール
6	0905-1420	PE シール(2 個入)
7	G4220-60016	バックアップリングを含めたシールホルダー
8	5062-2484	ガスケット、ウォッシュ用シール用(6 個入)
9	0905-1718	ウォッシュ用シール PE
10	G4220-63010	サポートリング (ウォッシュ用シール)
11		ポンプヘッドフェラル (ポンプヘッドをご注文ください)
12		ピストンハウジング (ポンプヘッドをご注文ください)
13	5067-5678	ピストン 1290 Infinity ポンプ(セラミック製)

10 メンテナンス用部品と器材 プライマリポンプヘッドの部品

プライマリポンプヘッド、シールウォッシュなし (バイナリポンプ)



メンテナンス用部品と器材 10 プライマリポンプヘッドの部品

品目	部品番号	説明
1	0515-1218	ネジ、M5、長さ 40 mm
2		ポンプチャンバハウジング (ポンプヘッドをご注文ください)
3	G4220-60022	インレットバルブ (プライマリポンプヘッド)
4	64220-60028	アウトレットバルブ 1290 Infinity クォータナリ ポンプ (プライマリポンプヘッド)
5	G4220-20020	1290 Infinity アウトレットバルブ用の内部ゴール ドシール
6	0905-1420	PE シール(2 個入)
7	G4220-60015	バックアップリングを含めたサポートリング
8		ポンプヘッドフェラル (ポンプヘッドをご注文ください)
9		ピストンハウジング (ポンプヘッドをご注文ください)
10	5067-5678	ピストン 1290 Infinity ポンプ (セラミック製)

10 メンテナンス用部品と器材 セカンダリポンプヘッドの部品

セカンダリホンノヘットの部品

セカンダリポンプヘッドの部品

セカンダリポンプヘッド、シールウォッシュ付き (バイナリポンプ)



図 22 セカンダリポンプヘッド、シールウォッシュ付 (バイナリポンプ)

メンテナンス用部品と器材 10 セカンダリポンプヘッドの部品

品目	部品番号	説明
1	0515-1218	ネジ、M5、長さ 40 mm
2	G4220-20003	ポンプヘッドネジ
3		ポンプヘッドフロントプレート (ポンプヘッドをご注文ください)
4	G4220-81013	熱交換器(セカンダリポンプヘッドのみ) チャンネル A
5	0515-5237	熱交換器ネジ
6	G4220-20001	スペーサフィッティング
7		ポンプチャンバハウジング (ポンプヘッドをご注文ください)
8	0905-1420	PE シール(2 個入)
9	G4220-60016	バックアップリングを含めたシールホルダー
10	5062-2484	ガスケット、ウォッシュ用シール用(6 個入)
11	0905-1718	ウォッシュ用シール PE
12	G4220-63010	サポートリング (ウォッシュ用シール)
13	(2x)	ポンプヘッドフェラル (ポンプヘッドをご注文ください)
14		ピストンハウジング (ポンプヘッドをご注文ください)
15	5067-5678	ピストン 1290 Infinity ポンプ (セラミック製)
16	G4280-60026	高圧フィルタアセンブリ (セカンダリポンプヘッド)
17	01018-22707	PTFE フリット(5 個)
18	5067-4728	シールキャップ

10 メンテナンス用部品と器材 セカンダリポンプヘッドの部品

セカンダリポンプヘッド、シールウォッシュなし (バイナリポンプ)



図 23 セカンダリポンプヘッド、シールウォッシュなし (バイナリポンプ)

メンテナンス用部品と器材 10 セカンダリポンプヘッドの部品

品目	部品番号	説明
1	0515-1218	ネジ、M5、長さ 40 mm
2	G4220-20003	ポンプヘッドネジ
3		ポンプヘッドフロントプレート (ポンプヘッドをご注文ください)
4	G4220-81013	熱交換器(セカンダリポンプヘッドのみ) チャンネル A
5	0515-5237	熱交換器ネジ
6	G4220-20001	スペーサフィッティング
7		ポンプチャンバハウジング (ポンプヘッドをご注文ください)
8	0905-1420	PE シール(2 個入)
9	G4220-60015	バックアップリングを含めたサポートリング
10	(2x)	ポンプヘッドフェラル (ポンプヘッドをご注文ください)
11		ピストンハウジング (ポンプヘッドをご注文ください)
12	5067-5678	ピストン 1290 Infinity ポンプ(セラミック製)
13	G4280-60026	高圧フィルタアセンブリ (セカンダリポンプヘッド)
14	01018-22707	PTFE フリット(5 個)
15	5067-4728	シールキャップ

- **10 メンテナンス用部品と器材** パージバルブ
- パージバルブ



図 24 パージバルブの部品

品目	部品番号	説明
	5067-4119	パージバルブヘッド
1	1535-4857	ステータネジ
2	5068-0004	パージバルブステータ
3	5068-0005	パージバルブのローターシール、1200 bar
4	1535-4045	ベアリングリング

メンテナンス用部品と器材 10 カバー部品

カバー部品



図 25 カバー部品

品目	部品番号	説明
1	5067-4613	キャビネットキット(左右の側面カバー、上部カ バー、チューブプラグ、下部カバー、リークシール)
2	5042-9949	チューブプラグ、プラスチック製
	5042-9972	チューブグロメット
3	5067-4612	前面カバー 1290 Infinity バイナリポンプ
4	5067-4634	バルブレールキット

10 メンテナンス用部品と器材 リーク部品

リーク部品



図 26	漏斗
------	----

部品番号	説明
5041-8388	漏斗

ヒューズ

部品番号		ĩ	兑明			
2110-1004	(2x)	ł	ニューズ	10	А	t

アクセサリキット

アクセサリキット

部品番号	説明
G4220-68705	アクセサリキット
5042-9974	チューブフレックス(1.5 m)
8710-0510	1/4 インチ × 5/16 インチのレンチ
8710-1924	両ロスパナ 14 mm
5023-0240	六角ドライバ、¼ インチ、スリット入り
8710-2392	六角レンチ4 mm15 cm 、T 字型ハンドル
9301-0411	シリンジ (プラスチック製)
9301-1337	シリンジアダプタ
5067-4657	SST キャピラリ 300 x0.17 mm、ポンプからオートサン プラ間
5067-4658	SST キャピラリ 450 x0.17 mm、ポンプからサーモス タット付きオートサンプラまで
0100-1710	チューブ接続用取り付けツール
G4220-67000	廃液チューブ
8710-1534	レンチ、4 mm 両端、オープンエンド
5065-9978	チューブ、内径1 mm 、外径3 mm 、シリコン、5 m、 ウォッシュ用シールオプション用
5181-1519	CAN ケーブル、Agilent モジュール間、1 m
01018-22707	PTFE フリット(5 個)
01018-23702	インサートツール
5042-9972	チューブグロメット

10 メンテナンス用部品と器材 その他

その他

品目	部品番号	説明
	5067-4699	1290 Infinity ポンプサービスキット
1	G4220-20012	トルクレンチ 2 - 25 Nm
2	G4220-20013	4 mm 六角ビット
3	G4220-20014	2.5 mm 六角ビット
4	G4220-20015	アダプタ ¼ インチ、正方形 / 六角形
	G4220-44000	保護用緩衝材
	5023-0285	1290 Infinity ポンプヘッド アライメント ツー ル(ピストン / ハンドル)用交換キット
	G4203-68708	HPLC システムツールキット

1290 Infinity ポンプサービスキット (5067-4699) には、ポンプヘッドア ライメントツールとアイテム 1 - 4 が含まれます。



メンテナンス用部品と器材 10 その他



1290 Infinity ポンプヘッド アライメント ツール (ピストン/ハ ンドル) 用交換キット 図 28



図 29 HPLC システムツールキット

その他



Agilent 1290 Infinity バイナリポンプ ユーザーマニュアル

11 ケーブルの識別

ケーブル概要 214 アナログケーブル 216 リモートケーブル 218 BCD ケーブル 222 CAN/LAN ケーブル 225 RS-232 ケーブルキット 226 Agilent 1200 モジュールからプリンタ 227

本章では、すべてのケーブルに関する情報を記載します。



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」218 ページ 図』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 +
F			
- TE			



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	高
	5,14	7 - 赤	レディ	高
	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 III/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	高
	14	7 - 赤	レディ	高
	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
	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)
A 🔘 1	白	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
s 0 15	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
1. Star	紫	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)

11 ケーブルの識別

RS-232 ケーブルキット

RS-232 ケーブルキット

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

ケーブルの識別 11

Agilent 1200 モジュールからプリンタ

Agilent 1200 モジュールからプリンタ

部品番号 説明

5181-1529 ケーブル「プリンタシリアルおよびパラレル」は SUB-D 9 ピンのメスであるのに対して、もう一方はセントロニクスコ ネクタ (ファームウェア更新には使えません)です。G1323 コントロールモジュール用です。

11 ケーブルの識別

Agilent 1200 モジュールからプリンタ



Agilent 1290 Infinity バイナリポンプ ユーザーマニュアル

12 ハードウェア情報

ファームウェアについて 230 電気的接続 233 モジュールの背面図 234 インタフェース 235 インタフェースの概要 239 8 ビットコンフィグレーション スイッチの設定 243 特別な設定 247 機器レイアウト 249 EMF (Early Maintenance Feedback) 250

この章では、ハードウェアと電子機器に関してポンプの詳細を説明 します。



ファームウェアについて

本装置のファームウェアは、次の 2 つの独立したセクションで構成されて います。

- ・ **レジデントシステム**と呼ばれる機器固有ではないセクション
- ・ メインシステムと呼ばれる機器固有のセクション

レジデントシステム

ファームウェアのレジデントセクションは、すべての Agilent 1100/1200/1220/1260/1290 シリーズモジュールで同一です。次のような機 能があります。

- 全通信機能 (CAN、LAN、および RS-232C)
- メモリー管理
- 「メインシステム」のファームウェアを更新する機能

メインシステム

次のような機能があります。

- 全通信機能 (CAN、LAN、および RS-232C)
- メモリー管理
- ・「レジデントシステム」のファームウェアを更新する機能

この他にメインシステムが備えている機器機能は、次のような一般機能に 分類できます。

- APG リモートを経由した同期実行
- エラー処理
- 診断機能
- 次のモジュール特有の機能
 - ランプコントロール、フィルタ動作、
 - 生データ収集、吸光度への変換などの内部イベント。

ファームウェアの更新

ファームウェアの更新は、以下のユーザインターフェイスから行うことができます。

- ハードディスク上のローカルファイルを用いた PC とファームウェアの 更新ツール
- USB フラッシュディスクのファイルを用いたインスタントパイロット (G4208A)
- Agilent LabAdvisor ソフトウェア(B.01.03 以降)

ファイル名の付け方は、次の規則に従っています。

PPPP_RVVV_XX.dlb、ここで

PPPP は製品番号です。たとえば、G1315A/B DAD の 1315AB です。

R はファームウェアの改訂のことです。たとえば、G1315B の場合は A、G1315C DAD の場合は B です。

WW は、改訂番号です。たとえば、102 はリビジョン 1.02 です。

XXX はファームウェアのビルド番号です。

ファームウェアの更新の説明については、メンテナンスの章のファーム ウェアの置換のセクション、またはファームウェアの更新ツールのドキュメ ントを参照してください。

注記 メインシステムの更新は、レジデントシステムにおいてのみ可能です。レジデ ントシステムの更新は、メインシステムにおいてのみ可能です。

メインシステムとレジデントシステムは同じセットのものである必要がありま す。 ファームウェアについて



図 30 ファームウェア更新の仕組み

注記

一部のモジュールは、そのメインボードのバージョンや初期ファームウェアバージョンにより、ダウングレードに制限があります。たとえば、G1315C DADSLをファームウェアのリビジョン B.01.02 以前や A. xx. xx にダウングレードすることはできません。

モジュールの中には特定のコントロールソフトウェア環境での操作を可能にす るために復旧できるものがあります(G1314CからG1314Bなど)。この場合、 復旧後のタイプの機能セットは使用できますが、復旧前の機能セットは失われ ます。再度、復旧処理を行うと(G1314BからG1314Cなど)オリジナルの機 能セットが再び使用できるようになります。

これら具体的な情報のすべては、ファームウェアの更新ツールのドキュメントに記載されています。

ファームウェアの更新ツール、ファームウェア、ドキュメントは Agilent のウェブサイトから入手できます。

 http://www.chem.agilent.com/EN-US/SUPPORT/DOWNLOADS/FIRM-WARE/Pages/LC.aspx

電気的接続

- CAN バスは、高速データ転送機能を持つシリアルバスです。CAN バスの
 2 つのコネクタは内部モジュールのデータ転送および同期に使用されます。
- スタートや、ストップ、共通シャットダウン、プレランなどの機能を利用したい場合は、リモートコネクタを他の Agilent Technologies 製分析機器と組み合わせて使用してください。
- 適切なソフトウェアを使用すれば、RS-232C コネクタを使って、コン ピュータから RS-232C 接続を介してモジュールをコントロールすること ができます。このコネクタは、コンフィグレーション スイッチで有効に し、設定することができます。
- 電源ケーブルコネクタは、100 240 VAC ± 10 % の入力電圧、50 または 60 Hz の電源周波数に対応しています。最大消費電力はモジュールごとに異なります。パワー サプライは広範な対応機能を備えているので、モジュールには電圧スイッチがありません。また、パワー サプライには自動電子ヒューズが装備されているため、外部のヒューズは必要ありません。
- 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。

12 ハードウェア情報

電気的接続





電源コネクタ

図 31 バイナリポンプの背面

インタフェース

Agilent 1200 Infinity シリーズのモジュールは、次のインターフェイス を装備しています。

モジュール	CAN	LAN/BCD (オプ ション)	LAN (オン ボード)	RS-232	アナログ	APG リモート	特殊
ポンプ							
G1310B Iso Pump G1311B Quat Pump G1311C Quat Pump VL G1312B Bin Pump G1312C Bin Pump VL 1376A Cap Pump G2226A Nano Pump G5611A Bio-Inert $ク オ - タ + U ポン \mathcal{T}$	2	あり	なし	あり	1	あり	
G4220A/B Bin Pump G4204A Quat Pump	2	なし	あり	あり	なし	あり	CAN スレー ブ用 CAN DC 出力
G1361A Prep Pump	2	あり	なし	あり	なし	あり	CAN スレー ブ用 CAN DC 出力

表 9 Agilent 1200 Infinity シリーズインターフェイス

インタフェース

表 9 Agilent 1200 Infinity シリーズインターフェイス

モジュール	CAN	LAN/BCD (オプ ション)	LAN (オン ボード)	RS-232	アナ ログ	APG リモート	特殊
サンプラ							
G1329B ALS G2260A Prep ALS	2	あり	なし	あり	なし	あり	G1330B 用サーモス タット
G1364B FC-PS G1364C FC-AS G1364D FC-mS G1367E HiP ALS G1377A HiP micro ALS G2258A DL ALS G5664A Bio-Inert FC-AS G5667A Bio-Inert オートサンプラ	2	あり	なし	あり	なし	あり	G1330B 用サーモス タット CAN スレーブ用 CAN DC 出力
G4226A ALS	2	あり	なし	あり	なし	あり	
検出器							
G1314B VWD VL G1314C VWD VL+	2	あり	なし	あり	1	あり	
G1314E/F VWD	2	なし	あり	あり	1	あり	
G4212A/B DAD	2	なし	あり	あり	1	あり	
G1315C DAD VL+ G1365C MWD G1315D DAD VL G1365D MWD VL	2	なし	あり	あり	2	あり	
G1321B FLD G1362A RID	2	あり	なし	あり	1	あり	
G4280A ELSD	なし	なし	なし	あり	あり	あり	外部接点 自動ゼロ

ハードウェア情報 12 インタフェース

モジュール	ュール CAN LAN/BCD LAN (オプ (オ ション) ボー		LAN (オン ボード)	RS-232	アナ ログ	APG リモート	特殊	
その他								
G1170A Valve Drive	2	なし	なし	なし	なし	なし	オン K LAN (フェア A) ルウェ A (フェア A) B. 06. 40 C. 06. 40 G4212A 40 G4212A 40 G4212A 50 G4212A 50 に G4212A 50 に 50 た 50 に 50 に 50 に 50 に 50 た 50 に 50 こ 5 こ 50 こ 50 こ こ 5 こ 5 こ 5 こ 5 こ 5 こ 5 こ 5 こ 5 こ	
G1316A/C TCC	2	なし	なし	あり	なし	あり		
G1322A DEG	なし	なし	なし	なし	なし	あり	AUX	
G1379B DEG	なし	なし	なし	あり	なし	あり		
G4225A DEG	なし	なし	なし	あり	なし	あり		
G4227A フレックスキューブ	2	なし	なし	なし	なし	なし		
G4240A チップキューブ	2	あり	なし	あり	なし	あり	CAN スレー ブ用 CAN DC 出力 G1330A/B 用 サーモス タット(不 使用)	

表 9	Agilent	1200	Infinity	シリーズイ	ンターフェイン	ス
-----	---------	------	----------	-------	---------	---

インタフェース

注記 LAN 経由での制御には、検出器 (DAD/MWD/FLD/VWD/RID) が望ましいアクセス ポイントとなります。モジュール間通信は、CAN を介して行います。

- CAN コネクタ (他のモジュールへのインターフェイス)
- LAN コネクタ (コントロールソフトウェアへのインターフェイス)
- RS-232C (コンピュータへのインターフェイス)
- リモートコネクタ(他のアジレント製品へのインターフェイス)
- アナログ出力コネクタ(シグナル出力用)

インタフェースの概要

CAN

CAN は、モジュール間通信インターフェイスです。これは、高速データ通信とリアルタイム要求をサポートする2線式シリアルバスシステムです。

LAN

注記

これらのモジュールには、LAN カード用インターフェイススロット (Agilent G1369B/C LAN インターフェイス)またはオンボード LAN イン ターフェイス(検出器 G1315C/D DAD や G1365C/D MWD など)が装備され ています。このインターフェイスにより、PC で適切なコントロールソフト ウェアを使用して、モジュール/システムを制御できます。

 Agilent 検出器 (DAD/MWD/FLD/VWD/RID) を使用したシステムの場合、LAN は DAD/MWD/FLD/VWD/RID に接続してください (データ負荷が高いため)。
 Agilent 検出器がシステムに含まれていない場合、ポンプまたはオートサンプ ラに LAN インターフェイスを取り付けてください。

RS-232C(シリアル)

RS-232C コネクタは、適切なソフトウェアを使用して、コンピュータから RS-232C 接続を介してモジュールをコントロールする場合に使用します。 このコネクタは、モジュールの背面にあるコンフィグレーションスイッチ モジュールで設定することができます。RS-232C の通信設定 を参照してく ださい。

注記 オンボード LAN を備えたメインボードで設定できるコンフィグレーションは ありません。これらは、あらかじめ以下のように設定されています。

- ボーレート 19200
- パリティなし 8 データビット
- スタートビット1つとストップビット1つは常に使用します(選択不可)。

RS-232C は、9 ピン (オス) SUB-D タイプコネクタを持つ DCE (データ通 信装置)として設計されています。ピンは次のように定義されています。

	表 1() RS-232	C 接続表
--	------	----------	-------

ピン	方向	機能
1	入力	DCD
2	入力	RxD
3	出力	TxD
4	出力	DTR
5		グランド
6	入力	DSR
7	出力	RTS
8	入力	CTS
9	入力	RI



RS-232 ケーブル 図 32

アナログシグナル出力

アナログシグナルは、記録用デバイスにも分配できます。詳細は、モ ジュールのメインボードの説明を参照してください。

APG リモート

他のアジレント製分析機器に一般的なシャットダウンや準備などの機能を 利用する場合、APG リモートコネクタを使用します。

リモートコントロールによって、シングル機器またはシステム間を簡単に 接続し、簡単なカップリング条件で、各機器を統合した分析が実行できま す。

リモートインターフェイスには、D コネクタを使用します。本モジュール は、入力 / 出力用 (ワイアード OR) リモートコネクタを 1 個装備してい ます。

各分析システム内での安全性を確保するために、1本はいずれかのモジュールで重大な問題が検出された場合に行うシステムの重要部分の[シャットダウン]専用になっています。すべての関連するモジュールがオンになっている(または正しく電源投入されている)ことを検出するため に、ラインの1本は接続されたすべてのモジュールの[電源オン]を要約するために定義されます。次の分析の準備を指示する[レディ]シグナル、その後、それぞれのラインで引き起こされる分析の[スタート]シグ ナルと[ストップ]シグナル(オプション)によって分析のコントロールを続けることができます。さらに、[プリペア]と[スタートリクエスト]も使用できます。シグナルレベルは次のように定義されています。

- ・ 標準 TTL レベル (0 V ロジック真、+ 5.0 V が偽)
- ファン出力は 10
- 入力負荷は 5.0 V に対して 2.2 kOhm
- ・ 出力はオープンコレクタ型、入力 / 出力 (ワイアード OR)

注記

一般的な TTL 回路はすべて、5 V パワーサプライで動作します。TTL シグナ ルは、0 V ~ 0.8 V の場合「低」または L、2.0 V ~ 5.0 V の場合「高」ま たは H と定義されます (それぞれ、アース端子に対して)。

表 11	リモー	・トシグナル	<i>、</i> ディスト	リビュ	ーション
------	-----	--------	---------------	-----	------

ピン	シグナル	説明
1	DGND	デジタルグランド
2	PREPARE	(L) 分析を準備するように要求します(キャリブ レーション、検出器ランプ点灯等)。受信側は、分析 前の動作を実行する任意のモジュールです。
3	START	(L) 測定 / タイムテーブルを開始するように要求しま す。受信側は、分析時間をコントロールできる任意 のモジュールです。
4	SHUT DOWN	(L)システムの重大な問題の発生を出力します(リークの発生時に ポンプを停止するなど)。ポンプ 停止)。受信側は、安全リスク軽減機能を持つ任意の モジュールです。
5		未使用
6	POWER ON	(H)システムに接続されたすべてのモジュールが ON になっていることを出力します。受信側は、他のモ ジュールの動作に依存する任意のモジュールです。
7	READY	(H)システムが次の分析の準備を完了していることを出力します。受信側は、任意のシーケンスコントローラです。
8	STOP	 (L)できるだけ早くシステムをレディ状態にするよう に要求します(測定の停止、注入の中断または終了)。受信側は、分析時間をコントロールできる任意の モジュールです。
9	START REQUEST	(L) インジェクションサイクルを開始するように要求 します (任意のモジュールでスタートキーが押され た場合等)。受信側はオートサンプラです。

特殊インターフェイス

モジュールは、例えば外部バルブなどの CAN スレイブとして動作するモジュールを共に使用することを目的とした DC-Out (24 VDC) 電力ライン を含みます。ラインは、0.5 A(1.7 A as of August 2011)の出力に制限され、自動リセットします。

8 ビットコンフィグレーション スイッチの設定

8 ビットコンフィグレーション スイッチの設定

8 ビットコンフィグレーションスイッチは、モジュール背面にあります。 このスイッチを使用して、LAN、シリアル通信プロトコル、機器固有の初期 化手順を指定するコンフィグレーションパラメータを設定できます。 オンボード LAN を搭載したすべてのモジュールの場合(G1315/65C/D、 G1314D/E/F、G4212A/B、G4220A/B など):

- デフォルトはすべてのスイッチがダウン(最適な設定)です。
 - LAN 用の Bootp モード
 - RS-232 用ボーレート 19200、パリティなし 8 データビット
- 特殊な LAN モードの場合、必要に応じて、スイッチ3~8 を設定する 必要があります。
- BOOT/テストモードの場合、スイッチ1と2をアップすることに加え、 必要なモードに設定する必要があります。

通常の操作ではデフォルト(最適)設定を使用します。

注記

8 ビットコンフィグレーション スイッチの設定



図 33 コンフィグレーションスイッチの位置 (例は G4212A DAD)

8 ビットコンフィグレーション スイッチの設定

注記

LAN コンフィグレーションを行うには、SW1 および SW2 を OFF に設定する必要があります。LAN 設定/コンフィグレーションの詳細は、「LAN コンフィグレーション」の章を参照してください。

表 12 8 ビット コンフィグレーションスイッチ (オンボード LAN あり)

	モード		機能					
	SW 1	SW 2	SW 3	SW 4	SW 5	SW 6	SW 7	SW 8
LAN	0	0	リンク コン	フィグレー	ーション	初期	化モードi	選択
自動ネゴシ	エーショ	ン ン	0	Х	х	Х	Х	Х
10 MBit、ハーフ	'デュプレ	ックス	1	0	0	Х	Х	Х
10 MBit、フル	デュプレジ	ックス	1	0	1	Х	Х	х
100 MBit、ハーフ	フデュプレ	(ックス)	1	1	0	Х	х	
100 MBit、フル	デュプレ	ックス	1	1	1	X X		Х
Вос	otp		Х	х	х	0 0 0		0
Bootp お	よび保存		Х	х	х	0 0		1
保存されたパラ	ラメータを	使用	Х	Х	х	0 1		0
DH	CP		Х	х	х	1 0		0
デフォル	トを使用		Х	х	х	0 1 1		1
テスト	1	1	システム					NVRAM
Boot レジデ	ントシスラ	FД	1					х
デフォルトラ (コールド	データに戻 スタート	[す)	Х	Х	Х			1

8 ビットコンフィグレーション スイッチの設定

凡例:

0 (スイッチダウン)、1 (スイッチアップ)、x (任意の位置)

注記	TEST モードを選択した場合、LAN 設定は、「自動ネゴシエーション」および
	「保存されたパラメータを使用」です。

注記	「Boot レジデントシステム」および「デフォルトデータに戻す (コールドス タート)」の詳細についてけ 『「焼別な設定」 247 ページ 図』を参照してくだ
	クロトリーの評判については、「特別な設定」241、シン区」を参照してくたさい。

8 ビットコンフィグレーション スイッチの設定

特別な設定

固有の処理には特別な設定が必要です(通常はサービス事例で)。

注記

表は、オンボード LAN を装備した場合、装備していない場合の両方のモジュールの設定を示しています。それぞれを LAN および LAN なしと識別して表示します。

Boot - レジデント

ファームウェアローディングエラー(メインファームウェア部分)が発生 した場合、ファームウェア更新手順でこのモードが必要となることがあり ます。

以下のスイッチ設定を使用し、機器の電源を再び入れると、機器ファーム ウェアはレジデントモードを維持します。この場合、モジュールとして操 作することはできません。オペレーティングシステムの基本機能(通信な ど)のみが使用できます。このモードでは、メインファームウェアを読み 込むことができます(更新ユーティリティを使用)。

表 13 Boot レジデント設定 (オンボード LAN)

モード選択	SW1	SW2	SW3	SW4	SW5	SW6	SW7	SW8
テスト /BOOT	1	1	1	0	0	0	0	0

8 ビットコンフィグレーション スイッチの設定

強制コールドスタート

強制コールドスタートを使用して、モジュールをデフォルトパラメータ設 定の定義済みモードにできます。

注意 データ損失 強制コールドスタートは、不揮発性メモリに保存されたメソッドとデータ をすべて消去します。ただし、キャリブレーション設定と、診断および修 理ログブックだけは消去されずに保存されます。

→ 強制コールドスタートを実行する前に、メソッドおよびデータを保存してください。

スイッチを以下のように設定して、機器の電源を入れ直すと、強制コール ドスタートが完了します。

表 14 強制コールドスタート設定 (オンボード LAN)

モード選択	SW1	SW2	SW3	SW4	SW5	SW6	SW7	SW8
テスト /B00T	1	1	0	0	0	0	0	1

機器レイアウト

モジュールの工業デザインには、いくつかの革新的な特徴が含まれていま す。これは、電子装置と機械的アセンブリのパッケージングに関するアジ レントの E-PAC コンセプトに基づいています。このコンセプトの基本は、 発泡プラスチックスペーサの発泡ポリプロピレン(EPP)層を使用して、そ の中にモジュールのメカニカルボードおよびエレクトロニックボードコン ポーネントを納めることです。このパックが金属製内部キャビネットに組 み込まれ、さらにプラスチック外装キャビネットで覆われます。このパッ ケージ技術の利点として、以下のような点があります。

- 固定ネジ、ボルト、またはワイヤーを実際になくすことにより、コン ポーネント数が減り、取り付け/取り外しを素早く行うことができる。
- 冷却エアーが必要な位置に正確に導入されるように、プラスチック層内 にエアチャネルが成形されている。
- このプラスチック層は、物理的なショックから、電子部分と機械部分を 保護する。
- 金属製内部キャビネットによって、内部電子回路ボードを電磁妨害から 遮蔽し、機器自体からの無線周波放出を減少または排除する。

EMF (Early Maintenance Feedback)

EMF (Early Maintenance Feedback)

本機器のメンテナンスとして、機械的摩耗または応力にさらされる流路内 の部品を交換する必要があります。理想的には、部品を交換する頻度は、 あらかじめ決めた間隔ではなく、モジュールの使用頻度と分析条件に基づ いて決定します。Early Maintenance Feedback (Early Maintenance Feedback)機能は、機器内の各部品の使用状態をモニタリングし、ユー ザー設定可能なリミットを超えた時点でユーザーにフィードバックする機 能です。この機能は、ユーザーインターフェイスの表示によって、メンテ ナンス作業が必要な時期であることを知らせます。

EMF カウンタ

EMF カウンタ は、使用のたびに増分されます。カウンタの上限値を指定し ておき、その限度を超えた時点でユーザーインターフェイスにフィード バックすることができます。一部のカウンタは、必要なメンテナンス手順 の終了後にゼロにリセットできます。

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。 Agilent 1290 Infinity バイナリポンプ ユーザーマニュアル

LAN コンフィグレーション



13

最初の必要事項 252 TCP/IP パラメータコンフィグレーション 254 設定スイッチ 255 初期化モード選択 257 動的ホストコンフィグレーションプロトコル(DHCP) 262 一般情報 (DHCP) 262 セットアップ (DHCP) 264 リンクコンフィグレーション選択 266 Bootp を使用した自動コンフィグレーション 267 Agilent BootP サービスについて 267 BootP サービスの使用 268 状況: LAN 通信を確立できない場合 268 BootP サービスのインストール 269 MAC アドレスを決定する 2 つの方法 271 Agilent BootP サービスによる IP アドレスの割り当 T 273 Agilent BootP サービスによる機器の IP アドレスの変 更 276マニュアルコンフィグレーション 278 Telnet の使用 279 インスタントパイロット (G4208A) の使用 283 PC およびユーザーインターフェイス ソフトウェアの設 定 284 ローカル コンフィグレーションのための PC セットアッ プ 284 ユーザーインターフェイス ソフトウェアの設定 285

この章では、検出器の Agilent ChemStation PC への接続について 説明します。



13 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

モジュールにはオンボードの LAN 通信インタフェースが搭載されています。

 今後参照するために、MAC (Media Access Control) アドレスを書き留め てください。LAN インターフェイスの 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。別のネットワークデバイスが同 じハードウェアアドレスを持つことはありません。モジュールの背面の ラベルに MAC アドレスの表示があります(『253 ページ 図 35』を参照)。



ポンプメインボードの部品番号 リビジョンコード、ベンダー、 製造年と週 MAC アドレス 生産国

図 34 MAC ラベル
最初の必要事項

- **2** 機器の LAN インタフェース(『253 ページ 図 35』を参照)を
 - 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(ポイントツーポイント接続)、あるいは
 - ・標準 LAN ケーブルを使用してハブまたはスイッチに接続します。



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)」264ページ図』を参照)。
- Telnet を使用して、手動でパラメータを設定する。
- インスタントパイロット(G4208A)を使用してパラメータをマニュアル 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。『257 ページ表 16』を参照してください。

LAN コンフィグレーション 13 設定スイッチ

設定スイッチ



設定スイッチはモジュール背面にあります。

図 36 設定スイッチの位置

上記のように、モジュールはすべてのスイッチが OFF に設定されて出荷さ れます。

注記

LAN 設定を行うには、SW1 および SW2 を OFF に設定する必要があります。

設定スイッチ

表 15 出荷時の初期設定

初期化(「Init」)モード	Bootp、すべてのスイッチがダウン。詳細につ いては、『「初期化モード選択」257 ページ 図』 を参照してください。
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、詳細については、 『「リンクコンフィグレーション選択」266 ペー ジ 図』を参照してください。

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 16 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。



Agilent 1290 Infinity バイナリポンプ ユーザーマニュアル

LAN コンフィグレーション 13 初期化モード選択 **13** LAN コンフィグレーション 初期化モード選択

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。



注記

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。

初期化モード選択

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。



初期化モード選択

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。『260 ページ 表 17』を参照してください。



図 40 デフォルトを使用(原則)

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。

表 17 デフォルトパラメータを使用

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

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。

初期化モード選択

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 41 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)

- **必要なソフトウェ** 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 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]	Reload
DAI	D : DE64260019 (G1315D)	
Main Revision	B.06.41 [0002]	
Resident Revison	B.06.40 [0007]	Print
On-time	3d 01:33h	<u> </u>
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	
	T	Exit
Information on each m	odule.	10:08

図 42 インスタントパイロットの 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 18 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 19 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)」262ページ 図』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 20 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(『257 ページ 表 16』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 Settings	
BootP Tab File:	
C:\Documents and Setti	ngs\All Users\Application Data\Agilent\BootP\TabFile
Create Tab File	Edit BootP Addresses
Logging	
Do you want to log	bootP requests?
BootP Log File:	
C:\Documents and Se	ttings\All Users\Application Data\Agilent\BootP\LogFile
- Default Settings	
Subnet mask:	0.0.0
Calaman	
Gateway:	0.0.0
	OK Cancel Help

図 43 [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 44 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」276ページ 図』または『「Agilent BootP サービスの再起動」277ページ 図』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 インターフェイス カードラベルを使った MAC アドレスの直 接入力

- 1 機器の電源を切ります。
- 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。
 『252 ページ 図 34』と 『253 ページ 図 35』を参照。
- 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	

図 45 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」276ページ 図』または『「Agilent BootP サービスの再起動」277ページ 図』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 46 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。

LAN コンフィグレーション 13 Bootp を使用した自動コンフィグレーション

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

dit BootP Addresse	25				×
Hardware Address 006000111999	Host Name AgilentLC1	IP Address 10.1.1.101	Comment Agilent LC1 right	Subnet Mask 255.255.255.0	Gateway 0.0.0.0
005000222888	AgilentLC2	101.1.1.102	Agilent LC2 left	255.255.255.0	0.0.0.0
	_			1	
Add	<u>M</u> od	fy <u>D</u> el	etek	se <u>H</u>	elp



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(『276ページ 図 46』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。

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。





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 49 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(『「設定スイッチ」 255ページ 図』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 50 モジュールとの接続が確立されます。

マニュアルコンフィグレーション

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 >?	G1315C PP00000024	
command syntax	description	
? / sm <x.x.x.x> gw <x.x.x.x> exit ></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 51 Telnet コマンド

表 21 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\	system32\cmd.exe - telnet 134.40.27.95
>/ LAN Status Pag	re
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 52 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\	system32\cmd.exe - telnet 134.40.27
>ip 134.40.27.	.99
LAN Status Pag	ge
MAC Address	: 0030D30A0838
Init Mode	: Using Stored
TCP/IP Propert - active -	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
凶 53 Ielne	et - 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 **13 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

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 54 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。

インスタントパイロット(G4208A)の使用

モジュールをネットワークに接続する前に TCP/IP パラメータを設定する には、インスタントパイロット (G4208A) を使用できます。

- 1 初期画面から、[More] ボタンを押します。
- **2** [Configure] を選択します。
- 3 DAD ボタンを押してください。
- 4 LAN 設定までスクロールダウンします。

Configure - DAD		
		Edit
Setting	Value	
Symbolic Name	<not set=""></not>	<u>⊳₩₩</u>
Temperature Control	ON	Bal.
UV-Lamp Tag	Use UV-lamp anyway	
Cell Tag	Use cell anyway	
Analog Out 1	0V - 1V output range	
Analog Out 2	0V - 1V output range	
UV lamp	Stays off at power on	
VIS lamp	Stays off at power on	
LAN IP	134.40.27.95	
LAN Subnet Mask	255.255.248.0	
LAN Def. Gateway	134.40.24.1	Evit
	_	
		13:26
System Contro	oller DAD	
		_ \ _

- 図 55 インスタントパイロット LAN コンフィグレーション
- 5 [編集] ボタン(編集モード以外の場合のみ表示)をクリックし、必要な 変更を行い、[完了] ボタンをクリックします。
- 6 [終了]をクリックして画面を終了します。

PC およびユーザーインターフェイス ソフトウェアの設定

PC およびユーザーインターフェイス ソフトウェアの設定

ローカル コンフィグレーションのための PC セットアップ

この手順では、ローカルコンフィグレーションのモジュールのデフォルト パラメータと一致させるためにご使用の PC の TCP/IP 設定を変更する方法 について説明します(『「初期化モード選択」257 ページ 図』を参照)。

固定 IP アドレス		
Local Area Connection Properties	Internet Protocol (TCP/IP) Properties
General	General	
Connect using: IIII Xircom CreditCard Ethernet 10/100 + Modem 56	You can get IP se this capability. Oth the appropriate IP	tings assigned automatically if your network supports erwise, you need to ask your network administrator for settings.
Configure Components checked are used by this connection:	Use the folk Use the folk IP address Subnet mask: Default gatewa	address automatically wing IP address 192 . 168 . 254 . 1 255 . 255 . 248 . 0 y:
Install Uninstall Properties Description Transmission Control Protocol/Internet Protocol. The default wide area network protocol that provides communication across diverse interconnected networks. ✓ Show icon in taskbar when connected OK Cancel	C Ogtain DNS C Use the folk Preferred DNS Alternate DNS	Internet Protocol (TCP/IP) Properties ? General ? You can get IP settings assigned automatically if your network supports this capability. Otherwise, you need to ask your network administrator for the appropriate IP settings. Obtain an IP address automatically Use the following if address: IP address Subhert mask
DHCP 経由の自動 IP アドレス (特別なネットワーク IT 設定が必要)		Defadit geteway:

図 56 PC の TCP/IP 設定の変更

PC およびユーザーインターフェイス ソフトウェアの設定

ユーザーインターフェイス ソフトウェアの設定

付属のユーザーインターフェイスソフトウェアの設定ガイドに従って、 ユーザーインターフェイスソフトウェアをインストールします。

13 LAN コンフィグレーション PC およびユーザーインターフェイス ソフトウェアの設定



Agilent 1290 Infinity バイナリポンプ ユーザーマニュアル

14 付録

安全に関する一般的な情報 288 廃電気電子機器(WEEE)指令(2002-96-EC) 291 無線妨害 292 騒音レベル 293 アジレントのウェブサイト 294

この章では、安全性、法律、ウェブに関する追加情報を記載してい ます。



14 付録

安全に関する一般的な情報

安全に関する一般的な情報

安全記号

表	22	安全記号
1	~~	シェルワ

記号	説明
	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。
\$	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。
Â	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

警告は、

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。
付録 14

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。

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。

Agilent 1290 Infinity バイナリポンプ ユーザーマニュアル

14 付録

安全に関する一般的な情報

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。

付録 14

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、廃液電気および電子機器 (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。

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、地元の Agilent 営業所にお問い合わせ頂くか、 詳細については Agilent のホームページ (www.agilent.com) を参照してくだ さい。

無線妨害

安全規準または EMC 規格に適合した方法で装置を正しく動作させるため に、Agilent Technologies 製以外のケーブルは使用しないでください。

テストと測定

選別していない機器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したりする場合、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。

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 Lp 70dB (A) 未満
- 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

アジレントのウェブサイト

アジレントのウェブサイト

製品およびサービスの最新情報を知るには、以下の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

Agilent 1290 Infinity バイナリポンプ ユーザーマニュアル

索引

Α

Agilent Lab Advisor ソフト ウェア 82 Agilent Lab Advisor 82 Agilent 診断用ソフトウェ ア 82 Agilent ユーザー インターフェイ ス ソフトウェアの設 定 284 APG リモート 241

В

BCD 222 ケーブル Bootp サービス 停止 276 インストール 269 再起動 277 設定 277 257 Bootp および保存 **258** 自動コンフィグレーショ ン 267 初期化モード 257 デフォルトの使用 260 保存されたパラメータの使 用 259

С

CAN 通信消失 89

CAN 239 ケーブル 225

D

DHCP セットアップ 264 一般情報 262

E

EMF early maintenance feedback 250

J

Jet Weaver 146

L

Lab Advisor ソフトウェ ア テスト 120 LAN Bootp および保存 258 Bootp を使用した自動コン フィグレーション 267 Bootp 257 設定スイッチ 255 C およびユーザー イン ターフェイス ソフトウェ アの設定 284 TCP/IP パラメータコン フィグレーション 254

telnet を使用したマニュ アルコンフィグレーショ ン 279 最初の必要事項 252 ケーブル 225 コンフィグレーショ 251 ン 初期化モード選択 257 デフォルトの使用 260 保存されたパラメータの使 用 259 マニュアルコンフィグレー ション 278 リンクコンフィグレーショ ン選択 266

М

MAC アドレス 確認 273 MAC アドレス 252

Ρ

PC およびユーザーインター フェイス ソフトウェアの設 定 284

R

RS-232C 239 ケーブル 226

索引

Т

TCP/IP パラメータコンフィグ レーション 254 telnet コンフィグレーショ ン 279

あ

アウトレットバルブ 141 アジレント インターネット上 294 アナログ ケーブル 216 アナログシグナル 240 圧力が下限値より下回りまし t-97 圧力が上限圧カリミットを超 えました 96 圧力センサー 交換 137 圧力センサのキャリブレー ションが誤っているか、せん さがありません 116 安全クラス | 289 安全 規格 19 一般的な情報 289 記号 288

こ

インターネット 294 インターフェイス 235 一般エラーメッセージ 86

え

エラーメッセージ 85

lost CAN partner 89 圧力が下限値より下回りま した 97 圧力が上限圧力リミットを 招えました 96 圧力センサのキャリブレー ションが誤っているか、せ んさがありません 116 オンにしようとしても ウォッシュ用シールポンプ がありません 117 カバー違反 95 カバーなしで起動 94. 94 シャットダウン 87 セカンダリポンプドライブ の電流が高過ぎま す 113 送液のアンダーラ ン 107 タイムアウト 86 デガッサの圧カリミット違 98 反 ドライブエンコーダ LED の電力が不十分 110 ドライブエンコーダエ ラー 111 ドライブエンコーダが機能 しなくなりました 104 ドライブタイムアウ ト 105 システム圧力テス 123 ト ドライブの電流が少なすぎ ます 104 ドライブの電流が高すぎま 105 す ドライブポジションリミッ 110 ト

パージバルブ故障 108 パージバルブのタグの書き 込み失敗 111 パージバルブのタグの読み 取り失敗 109 廃液カウンタの制限を超え ました 100 バイナリポンプが解析中に チャットダウン 102 バイナリポンプのデガッサ のターゲット圧力に達して いません 98 ファン動作不良 93 不明なパージバルブ タイ プ 114 プライマリポンプドライブ の電流が高過ぎま す 112 補正センサーオープ 92 ン 補正センサーショー ト 92 ポンプエンコーダの書き込 みが失敗しました 103 ポンプエンコーダの読み取 りが失敗しました 102 ポンプドライブが詰まった かまたはエンコーダが機能 しなくなりました 103 ポンプドライブエンコーダ のエラー 114 ポンプドライブエンコーダ の不具合 108 ポンプドライブエンコーダ のロールオーバー 109 ポンプドライブストップが 見つかりません 116 ポンプドライブストローク がブロックされまし た 115

ポンプヘッドのリークテス F 126 ポンプドライブのエ - 115 ポンプドライブの過電 流 106 ポンプのエラー 96 メインボードとポンプドラ イブエンコーダの接続不具 合 107 溶媒カウンタがリミットを 招えました 99 溶媒切り替えバルブ(SSV) の過電流 106 リークセンサーオープ 91 ン リークセンサーショー F 90 リーク 94 リモートタイムアウ ト 88 流速の制限を超えまし t-101

お

オンにしようとしてもウォッ シュ用シールポンプがありま せん 117 温度センサー 94

ታ

カラム外ボリューム 68 開梱 24

き

キャピラリ 134 機器レイアウト 249 凝縮 18

< クリーニング 133 け ケーブル BCD 222. 215 225. 215 CAN 225. 215 LAN RS-232 226. 215 アナログ 216. 214 214 概要 リモート 218. 214 C

コンフィグレーション 1 スタック 26 2 スタック 29 2 スタック前面 29 2 スタック背面 30 低流速 42 梱包チェックリスト 25 梱包の 傷み 24

さ

最小流速 42 最適化 67 高分解能達成 71 スタック コンフィグレー ション 26

し

システムの設定とインストー ル スタック コンフィグレー ションの最適化 26 シャットダウン 87

シールウォッシュオプショ ン 取り付け 40 シールウォッシュポンプ 交換 149 シャットオフバルブパネ ル 135 交換 シャットオフバルブ 交換 135 システム圧力テスト エラーメッセージ 123 設置について 15 溶媒キャビネット 38 仕様 15 湿度 19 周波数範囲 19 修理 ファームウェアの交 換 185 使用温度 19 使用高度 19 使用周囲温度 19 消費電力 19 仕様 物理的 19 初期化モード選択 257 診断用ソフトウェア 82 重量 19 真空脱気ユニット 42

す

ステータスインジケー タ 80 寸法 19

索引

せ

セカンダリポンプドライブの 電流が高過ぎます 113 設置スペース 18 設置要件 電源コード 17 設定スイッチ 255 設置 設置スペース 18 電源について 16

そ

送液のアンダーラン 107 た タイムアウト 86

τ

テスト 120 テスト機能 78 ディレイボリューム 68 説明 68 デガッサの圧カリミット違 反 98 電源周波数 19 電圧範囲 19 電気的接続 の説明 233 電源インジケーター 79 電源ケーブル 17 電源について 16

لح

トラブルシューティング エラーメッセージ 85, 78

ステータスインジケー ター 78 ドライブエンコーダ LED の電 力が不十分 110 ドライブエンコーダエ ラー 111 ドライブエンコーダが機能し なくなりました 104 ドライブタイムアウ ト 105 ドライブの電流が少なすぎま 104 す ドライブの電流が高すぎま す 105 ドライブポジションリミッ 110 ト 特殊インターフェイ 242 ス 特別な設定 強制コールドスター 248 ト Boot- レジデント 247 到着時不良 24

に

入力電圧 19

は

バイナリポンプが解析中に シャットダウン 102 バイナリポンプのデガッサの ターゲット圧力に達していま せん 98 バッファアプリケーショ ン 42 パージバルブの故障 108 パージバルブのタグの書き込 み失敗 111 パージバルブのタグの読み取 り失敗 109 パフォーマンス 最適化 67 廃液カウンタの制限を超えま 100 した 配管 ポンプへ 36 取り付け シールウォッシュオプショ ン 40 8 ビットコンフィグレーショ ンスイッチ オンボード LAN 243

ふ

ファームウェア アップグレード / ダウング レード 185 更新ツール 231 更新 185. 231 説明 230 メインシステム 230 レジデントシステ ム 230 ファン動作不良 93 フィッティング 134 フィルタ 溶媒インレット 42 プライマリポンプドライブの 電流が高過ぎます 112 不明なパージバルブ タイ プ 114 物理的仕様 19 部品 破損 25 不足 25

分解能

最適化 71

ほ

ポンプエンコーダの書き込み が失敗しました 103 ポンプエンコーダの読み取り が失敗しました 102 ポンプドライブが詰まったか またはエンコーダが機能しな くなりました 103 ポンプドライブエンコーダの エラー 114 ポンプドライブエンコーダの 不具合 108 ポンプドライブエンコーダの ロールオーバー 109 ポンプドライブストップが見 つかりません 116 ポンプドライブストロークが ブロックされました 115 ポンプドライブのエ ラー 115 ポンプドライブの過電 流 106 ポンプのエラーメッセー ジ 96 ポンプ 概要 11 準備 42 特徴 10 ポンプヘッドのリークテス ト エラーメッセージ 126 ポンプ 配管 36 原理 12

19

保管温度

保管高度 19 保管周囲温度 19 補正センサーオープン 補正センサーショート

マ

マニュアルコンフィグレー ション LAN 278

め

メインボードとポンプドライ ブエンコーダの接続不具 合 107 メッセージ カバー違反 95 カバーなしで起動 94. 94 リモートタイムアウ ト 88 メンテナンス 概要 128 ファームウェアの交 185 換 フィードバック 250

ŧ

藻 42

Ю

ユーザー インターフェイ ス 81

よ

溶媒インレットフィル タ 42

溶媒カウンタがリミットをこ えました 99 溶媒キャビネット 42 溶媒切り替えバルブ(SSV)の 92 過電流 106

IJ

92

リークセンサーオープ ン 91 リークセンサーショー F 90 リーク 94 リモート ケーブル 218 流速の制限を超えまし た 101 流速 42 最小 リンクコンフィグレーション 選択 266

www.agilent.com

本書の内容

本書には、Agilent 1290 Infinity バイナリ ポンプ G4220A に関する技術的リファレンス 情報が記載されています。

- 概要と仕様
- 設置、
- ・ 使用と最適化、
- トラブルシューティングおよび診断
- ・ メンテナンス
- 部品の識別、
- ハードウェア情報、
- ・ 安全保護と関連情報。

© Agilent Technologies 2010-2012

Printed in Germany 07/2012



G4220-96006 Rev. B

