Agilent 1290 Infinity LC システム

システムマニュアルおよび クイックリファレンス





Agilent Technologies



© Agilent Technologies, Inc. 2009-2011, 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

G4220-96301

エディション

05/2012

Printed in Germany

Agilent Technologies Hewlett-Packard-Strasse 8 76337 Waldbronn

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

保証

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

技術ライセンス

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

安全に関する注意

注意

注意は、危険を表します。これは、正しく実行しなかったり、指示を順守しないと、製品の損害または重要なデータの損失にいたるおそれがある操作手順や行為に対するたくかけたで、指示された条満たされるまで、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。

本ガイドの内容

本書では、Agilent 1290 Infinity LC システムについて説明します。

1 高速液体クロマトグラフィの紹介

本章では、Agilent 1290 Infinity LC システムとその基盤となっている概 念について紹介します。

2 Agilent 1290 Infinity LC システムの製品説明

この章では、1290 Infinity LC システムの特徴について説明します。

3 Agilent 1290 Infinity LC システムの最適化

この章では、理論の適用方法と、LC システムの特徴を活かして最適分離を 達成する方法について説明します。

4 システムの設定とインストール

この章では、ソフトウェアのインストール、スタック構成、システム運用 の準備について説明します。

5 クイックスタートガイド

この章では、1290 Infinity LC システムを使ったデータ取込とデータ分析 について説明します。

6 付録

この章では、安全性、法律、ウェブ、およびメソッドの設定に関する追加 情報を記載しています。 目次

1 高速液体クロマトグラフィの紹介 7 液体クロマトグラフィで、より小さな粒子を使用することの理論的根 拁 8 サブ2マイクロン粒子で充填カラムの利点 14 摩擦加熱 18 2 Agilent 1290 Infinity LC システムの製品説明 21 Agilent 1290 Infinity LC システムの新機能 22 システムコンポーネント 25 3 Agilent 1290 Infinity LC システムの最適化 39 ディレイボリュームとエクストラカラムボリューム 40 最適ディレイボリュームの設定方法 42 注入量を増加させる方法 51 ハイスループットを達成する方法 53 高分解能を達成する方法 56 感度を向上させる方法 59 キャリーオーバを最少にする方法 67 カラムの詰まりを回避する方法 69 4 システムの設定とインストール 71 ソフトウェアのインストール 72 モジュールの設置 74 5 クイックスタートガイド 91 クイックスタートガイドについて 92

システムの準備 93
 Method and Run Control (メソッド & ランコントロール) ビュー内の
 データ取得 98
 データ分析 107

4

6 付録 113
 安全に関する情報 114
 溶媒情報 117
 アジレントのウェブサイト 118
 メソッド全体の編集を使ったメソッドの設定 119

目次

目次



Agilent 1290 Infinity LC システムマニュアルおよびクイック リファレンス

高速液体クロマトグラフィの紹介

液体クロマトグラフィで、より小さな粒子を使用することの理論的根拠
サブ2マイクロン粒子で充填カラムの利点
14
摩擦加熱
18

本章では、Agilent 1290 Infinity LC システムとその基盤となって いる概念について紹介します。



液体クロマトグラフィで、より小さな粒子を使用することの理論的根拠

液体クロマトグラフィで、より小さな粒子を使用することの理論的根拠

はじめに

アジレントは、2003 年に、1.8 µm の粒子を使用する、最初の商用多孔質 シリカカラムを開発しました。これらのカラムは、後に「サブ 2 マイクロ ン」カラムまたは STM カラムとして知られるようになった最初の一連のカ ラムでした。これらの充填剤は、2006 年に開発された Agilent 1200 シ リーズ Rapid Resolution LC システムでの使用を目的として、600 bar の 耐圧力を持つ ZORBAX RRHT カラムで導入されました。2009 年には、 1200 bar の圧力でも日常的に使用可能な RRHD カラムまで含むように範囲 が拡大されました。これは最大で 1200 bar の圧力と 5 ml/min の流量の 運用能力を持つ新製品の Agilent 1290 Infinity LC システムをサポート するためでした。

このようなサブ 2 マイクロン (1.8 µm) の粒子径を持つカラムは、主として次の 2 つの目的で使われます。

1 クロマトグラフィの時間短縮

サブ 2 マイクロンの粒子サイズを持つ短いカラムは、分解能を損なうこ となく流量を増加させることができるため、分析に必要な時間を著しく 短縮します。

2 より優れた分解能

サブ 2 マイクロンの粒子サイズを持つ長いカラムは、効率が上がるため、複雑なサンプルの分離で必要とされる、より優れた分解能を獲得できます。また、特に LC/MS の場合は、拡散が少なくなるため、対象化合物のピークおよびその結果の感度のゲインの劣化が減少します。

STM 粒子を持つカラムを使うと、溶媒を促進するために必要な圧力を急激 に増加させることができます。これは分離の時間を短縮するために流量を 増加できること、そしてカラムを長くすればするだけ、分解能を向上させ ることができるからです。STM カラムが認知されたことによって、UHPLC システムを開発することが自然な流れとなりました。これは、HPLC の初期 の頃から続いてきた 400 bar の標準よりも高い圧力を提供できる HPLC シ ステムです。UHP (超高速または超高圧) LC システムは、クロマトグラ フィに高速高性能の分離性能を持たせるために必要とされる、ディレイボ

液体クロマトグラフィで、より小さな粒子を使用することの理論的根拠

リュームの低減とデータ収集の高速化に効果があります。Agilent 1290 Infinity LC は、UHPLC の分野での偉大な進歩です。市場に存在した各種 の UHPLC が持っていた各種の性能範囲をすべて完全にカバーする最初のシ ステムです。

理論

HPLC の分離効率は、van Deemter 方程式で表現できます(『9ページ 図 1』)。これは、対象化合物をカラムに沿って下げていく際に、拡散を測定 するために使われてきた段高さモデルの拡張です。H は理論段相当の高さ (HETP と表現されることもあります)、d_p はカラム充填剤の粒子径、u₀ は 移動相のリニア速度、そして A、B、C は各種の拡散力に関連付けられる定 数です。A 項は、渦拡散またはカラムを通過する複数の流路に関連付けら れます。B 項は、カラム軸(縦方向)に沿った分子拡散に関連付けられま す。C 項は、移動相と固定相間の対象化合物の物質移動に関連付けられま す。分離効率が最大になるのは、H が最小値のときです。各項とそれが結 合された方程式の効果を『9ページ 図 1』に示します。そこでは、プレー ト高さをカラム内の線型流量に沿ってプロットしています。この種のプ ロットは Van Deemter 曲線として知られており、カラムの分離効率を最大



液体クロマトグラフィで、より小さな粒子を使用することの理論的根拠

『10ページ 図 2』の van Deemter プロットでは、粒子径を小さくすること が効率を上げる結果となることが示されています。通常使われる 3.5 µm と 5.0 µm の粒子径を 1.8 µm の粒子サイズに変更すると、性能は著しく 向上します。1.8 µm の粒子では、プレート高さが 1/2 ~ 1/3 ほど低くな り、それに比例して効率が向上します。それに従って、分解能を損なうこ となくカラムを短くすることが可能になり、それに応じて分析時間も 1/2 ~ 1/3 ほど短縮されます。効率が向上するのは、粒子径を小さくした(つ まり、A 項(渦拡散)を小さくした)ために、複数の流路が短縮されたこ とが大きな理由です。また、粒子径を小さくすれば、物質移動の時間も短 縮できるため、C 項も小さくなります。そして、流量が増加した(線の傾 斜が緩やかになった)ため、全体の効果としては、効率損失が大幅に減少 したことが観察できます。つまり、流量を増加させればさせるほど、効率 の低下を招くことなく、小さな粒子の分離を加速できることがわかります。



隙間の線速度 (µ - cm/秒)

図 2 各種の粒子径に対する Van Deemter 曲線

Agilent 1290 Infinity LC システムマニュアルおよびクイックリファレンス

液体クロマトグラフィで、より小さな粒子を使用することの理論的根拠

クロマトグラフィの分離性能は、粒子径、孔径、粒子形態、カラムの長さ と内径、溶媒速度、温度のような、HPLC カラムの物理特性に基づいて最適 化できます。また、分離熱力学に配慮し、溶質と固定相 / 移動相の特性(有機溶媒の割合、イオン強度、および pH)を操作することによって、最小 の保留時間と最大の選択性を達成できます。



図 3 HPLC の最適条件の選択

分解能は、次の3 つのパラメータを持つ関数で表現できます。

- カラム効率または理論段数 (N)
- 選択性 (a)
- 保持係数(k)

分解能方程式(『11ページ図4』)によれば、分解能に与える影響は選択 性が最大であることがわかります(『12ページ図5』)。つまり、適切な 固定相/移動相の特性と温度を選択することが、分離を成功させるために 最も重要であるということです。



液体クロマトグラフィで、より小さな粒子を使用することの理論的根拠



図 5 R に対する段数、分離係数、保留係数の効果

高速分離メソッドは、新規に開発したものか、既存の汎用メソッドを単純 に流用したものかにかかわらず、各種のカラム形式で利用可能な固定相の 化学的性質の広範囲な選択を可能にするため役に立ちます。

アジレントは、すでに 140 種類以上の ZORBAX 1.8 µm 高速分離ハイス ループット(RRHT) カラム(14 の選択肢、15 ~ 150 mm の長さ、2.1、 3.0、4.6 mm の内径)を提供済みでしたが、Agilent 1290 Infinity LC の 開発に伴って、STM の範囲が高速分離ハイ デフィニション(RRHD) 1200 bar カラムまで含むように拡張されました。これによって、選択性を 最大化するために、最適の固定相が選択できるようになりました。分離性 能、流量、分析時間は、カラムの適切な長さと内径を選択することで最適 化できるため、長い STM カラムが従来よりも操作しやすくなりました。

多くのラボでは、分離用の固定相、移動相、温度の最適の組み合わせを選 択するために、念入りなスクリーニングプロセスを実施しています。アジ レントは、1200 シリーズと 1290 Infinity LC システムの双方に対して、 メソッド開発ソリューションを提供しています。これらのソリューション を使うと、時間がかかる選択プロセスを完全に自動化できます。そのため、 メソッド開発とメソッド移管が簡単で信頼性のある作業に変化します。

液体クロマトグラフィで、より小さな粒子を使用することの理論的根拠

ZORBAX 1.8 µm RRHD および RRHT カラムは、3.5 および 5 µm の粒子を使用する ZORBAX カラムと同じ化学的性質を使用します。その結果、任意の ZORBAX 相に対して、5.0、3.5、1.8 µm の粒子によって、同等の選択性が 提供されるため、従来の LC、UHPLC および準備中の LC 間で、簡単、高速 かつ安全な双方向のメソッド移管が実現できます。

サブ2マイクロン粒子で充填カラムの利点

サブ2マイクロン粒子で充填カラムの利点

クロマトグラフィの時間短縮

分析時間が短縮されることで、いくつかの利点がもたらされます。ハイス ループットのラボでは、キャパシティが増加するため、少ない時間でより 多くのサンプルを分析できるようになります。少ない時間でより多くのサ ンプルを分析できるため、コストが下がります。たとえば、分析時間をサ ンプルにつき 20 min から 5 min に短縮すると、700 個のサンプルのコス トは 79 % 削減されます (『14ページ 表 1』)。

表	1	700	分析の時間とコス	トの節減

サイクルタイム	20 分のサイクル タイム	5 分のサイクル タイム
分析	700	700
1 分析あたりの概算コスト ¹	\$ 10.58	\$ 2.24
700 分析あたりの概算コス ト ¹	\$ 7400	\$ 1570
コスト節減	_	\$ 5830
時間 ²	10 日	2.5 日

¹ 溶媒 = \$ 27/I 分析、廃棄処分 = \$ 2/I 分析、人件費 = \$ 30/時間

² 24 時間 / 日

Agilent のコスト節減カルキュレータには、従来の HPLC を 1.8 µm の粒 子径カラムを使った UHPLC に切り換えた場合のコスト節減高を計算する機 能があります。このカルキュレータは、メソッド移管カルキュレータと同 様に、アジレント テクノロジの Web サイト www.chem.agilent.com から 入手できます。結果はグラフ形式と表形式で表示されます。

分析時間が短縮されるということは、回答を得るまでの時間が短縮される ということでもあります。これは、プロセスコントロールと短時間のリ リーステストのために重要です。1 つの薬品の単一バッチをリリースする

サブ2マイクロン粒子で充填カラムの利点

ための、すべてのシステム適合性、キャリブレーション、サンプル分析の 実施に従来は数時間かかっていましたが、これが1時間以内に完了するよ うになります。短時間で回答を得られるということは、化合物の確認と反 応制御のためにオープンアクセス LC/MS システムを使用する合成化学者に とっても重要です。分析時間の短縮は、メソッド開発プロセスを著しく加 速します。

Agilent Technologies Method Translator and Cost Savings Calculator									
Basic Mode	Advanced N	lode	Viscosity	Table	Cost Savings Calculator	ן			
Settings							Financial View		
General Settings						^	Samples per Year 76500		
Solvent Costs (USD/liter)	3	.1	•				Max. Samples per Instrument pe	r Year (LC): 6480	
Waste Costs (USD/liter)	1	.4	•				Max. Samples per Instrument pe	r Year (RRLC): 19440	
Labour Costs (USD per year pe	r operator) 5	0000	•				600	10	
Depreciation in Y (linear calculation	rears 7		•				400	8	
Add. Laboratory (USD per year)	Costs)	•				200	4	
Daily Operating H (instrument)	lours 1	6	12 1	8 24			0	2	
Weekly Operating (instrument)	Days 1	2 3	4 5	6 7			Total Annual Operational Costs in Thousands of US	D	Cost per Sample in USD
Yearly Operating (instrument)	Weeks 1	10	20 30	40 50		=	Conventional LC	Agilent 1200 Series Rap	oid Resolution LC (RRLC) System
Instrument Setting	15	Convent	tional LC	Agilent 1200) Series RRLC System				
Instrument Costs (USD per instrume	ent)	50000		65000	•			Conventional LC	Agilent 1200 Series RRLC System
Uptime		90	A	90	÷		Operator Costs	590277.78 USD	196759.26 USD 👔
Maintenance Cos	ts	4000		4000			Maintenance Costs	47222.22 USD	15740.74 USD 🥑
(USD per instrume Column Costs	ent per year)	500		500	•		Instruments Costs	84325.40 USD	36541.01 USD 🕑
(USD per column)		500	•	500	•		Consumable Costs	0.00 USD	0.00 USD
(per instrument)		1	•	1	÷		Solvent Costs	5928.75 USD	411.87 USD 2
Consumable Cos (USD per sample)	ts 🕜	0	-	0	÷		Waste Costs	2677.50 USD	186.01 USD 👔
Method Settings							Annual Lab. Costs	0.00 USD	0.00 USD
Runtime 🕜 (min)		25	•	8.5	•				
Flow (ml/min)		1		0.21	÷		Total Annual Operational Co	768681.65 USD	287888.89 USD 🕡
Column Lifetime		 Inject 	tions				Total Cost per Sample	10.05 USD	3.76 USD 👔
		Col.	Volumes	1000	÷	•	Savings		480792.76 USD 🥑
						Currency	Switch to "Laboratory View		
Print Res	et Pro	ofiles	Help	Disclair	mer				Version 2.0 Online Update

図 6 コスト節減カルキュレータ

サブ2マイクロン粒子で充填カラムの利点



図 7 粒子径、効率、分析時間の間の関係

サブ2マイクロン粒子で充填カラムの利点

より優れた分離能

長いカラムでは、充填される粒子が小さければ小さいほど、効率と分離能 が向上します。このことは、メタボロミクスやプロテオミクスの研究のた めの複雑なサンプルの分析で重要です。また、不純物プロファイリングの ようなアプリケーションでは、高度な分離能力が役に立ちます。体液内の 薬品の LC/MS 分析でも、イオン抑制による干渉の減少で、ピークキャパシ ティの増加という利点が享受されます。一般に、分離能力が高ければ高い ほど、分析結果の信頼性が向上します。



図 8 BSA のトリプシン消化物の分析に ZORBAX RRHT SB-C18 カラム (2.1 x 150 mm、1.8 µm)を使った場合は、700 超のピークキャパシ ティを達成可能

1 高速液体クロマトグラフィの紹介 摩擦加熱



カラム内の移動相の圧力と流量を増加させていくと、熱が発生します。その結果発生する温度勾配(放射状および縦方向)は、カラムの効率に影響 を与えます。

Power = F * p

ここで、F は流量、p は圧力です。

強力なカラム温度調節機能(たとえば、水槽の使用)は、強い放射状の温 度勾配を生成し、それがカラム効率の著しい損失を招きます。静止空気に よるカラム温度調節機能は放射状温度勾配を平滑化するので効率の損失は 少なくなりますが、カラム出口から生ずる高い温度には耐性を持つ必要が あります。上がった温度は、選択性に影響します。背圧が低ければ低いほ ど摩擦加熱による性能損失が少なくなるため、内径が 4.6 または 3 mm の サブ 2 マイクロン カラムであっても、内径が 2.1 mm のカラムよりは、 効率が上です。

内径 2.1 mm の STM カラムへのグラジエント メソッド伝達で、分離速度 を上げていく場合の例を、『19 ページ 図 9』に示します。2.1 mm カラムの 最初の分析は、流量 0.22 ml/分、生成された圧力は 380 bar 、温度設定 は 37 °C 、すべてのピークは 12.5 分 ごとに分離された状態で実行され ています (クロマトグラムは非表示)。流量は 0.66 ml/分 に引き上げら れ、グラジエント時間は 1/3 に下方修正され、生成される圧力は 1020 bar で、すべてのピークが 4.2 分 ごとに現われるように変化してい ます (既出の『19 ページ 図 9』)。分離性能は同じであるべきですが、実 際にはピーク 7 と 8、およびピーク 5 とメインピークの間にカラム内の 加熱の影響による分解能の劣化が見られ、これらの化合物に対する選択性 を変化させています。カラム内の加熱効果を相殺し、分離性能を回復する には、カラムサーモスタットを 5 °C だけ下げれば十分であることがわか ります (下の『19 ページ 図 9』)。圧力は 1070 bar まで上がっています が、これもカラム内温度が低かったことを示しています。

高速液体クロマトグラフィの紹介 1 **摩擦加熱**



図 9 摩擦加熱発生が選択性とカラム温度低下の効果に与える影響

要約すれば、サブ 2 マイクロン充填剤の使用は、効率の向上、分解能の向上、分離時間の短縮という利点をもたらすことになります。Agilent 1290 Infinity LC システムと RRHD カラムによって、利用可能な分離空間が拡大されたため、上述の利点を利用できる機会が増大しました。1290 Infinity LC の特徴を『「Agilent 1290 Infinity LC システムの製品説明」 21ページ 図』に示します。また『「Agilent 1290 Infinity LC システムの 最適化」39ページ 図』では、これらの理論の適用方法とこれらの特徴を 活かして最適分離を達成する方法について説明します。

Agilent 1290 Infinity LC システムマニュアルおよびクイックリファレンス

19

1 高速液体クロマトグラフィの紹介 摩擦加熱 Agilent 1290 Infinity LC システムマニュアルおよびクイック リファレンス



2 Agilent 1290 Infinity LC システ ムの製品説明

Agilent 1290 Infinity LC システムの新機能22システムコンポーネント25

この章では、1290 Infinity LC システムの特徴について説明します。



Agilent 1290 Infinity LC システムの新機能

Agilent 1290 Infinity LC システムは、既存および今後出現する任意の種類のカラムテクノロジを使って液体クロマトグラフィ分析を実施する際に最大限の柔軟性を提供することを目的に設計されました。1290 Infinity LC は広範囲の運用パラメータに対応しているため、どのベンダーが製造した HPLC または UHPLC 分析システムであっても、インポートするメソッド設定の複製が可能です。Agilent 1290 Infinity LC システムは、この目標を達成するためのいくつかの革新的な設計コンセプトを備えており、これらの一連の優れた設計に裏付けられた信頼性によって、Agilent HPLC システムが、最も成功した HPLC ラインとして位置付けられています。

本システムは以下の特徴を備えています。

- 従来型のクロマトグラフィまたは高速クロマトグラフィに対応するため、0.05 ml/min~5 ml/min の流量、内径 1~5 mm のすべての分析カラム、そして任意のカラム充填剤をサポートします。
- ・最大許容圧力は 1200 bar (>17400 psi)です。そのため、サブ 2 マイ クロン充填剤と広範囲の移動相粘性選択肢を使って、短いカラムの場合 は高速クロマトグラフィ、長いカラムの場合は高い分解能を獲得できま す。
- ・ 質量分析検出機能または紫外可視吸光度検出機能を備えた、最高速グラ ジエントに対応した超低ディレイボリューム。
- 他の HPLC または UHPLC システムから移管する任意のメソッドを実行する機能。
- クロマトグラフノイズと音響ノイズを最小にするための高度なポンプコントロール機能。これによって、信頼性のある結果と働きやすい研究環境を獲得できます。
- ポンプモジュールに統合されたデガッサと自動パージバルブ。
- 低減したディレイボリューム、低減したキャリーオーバ、固定ループ オートサンプラとして動作するオプションを備えた可変ボリュームオー トサンプラ。
- オートサンプラに対して、ニードルシートのバック フラッシングや固定ループ動作のような機能を追加できる新規のフレキシブルキューブモジュール。

Agilent 1290 Infinity LC システムの新機能

- 使い勝手が強化されたカラム コンパートメントと、最大で 1200 bar (17400 psi)の圧力に対応した統合バルブソリューション。
- ・ 感度が著しく向上したダイオードアレイ検出器と、光学流体導波路を備 えたカートリッジセルシステムによるベースラインの安定性。
- ・ 全スペクトル情報を伴なった、最大 160 Hz のデータ取込レート。
- 一連の ZORBAX RRHD サブ 2 マイクロン粒子径カラム。動作最大圧力は、 1200 bar。
- 1290 Infinity クォータナリポンプの自動バッファリングと添加剤混合 に対応した混合アシスト。

最も重要な進化は、システムで使用できる圧力の範囲と流量が変わったこ とです。動作範囲は機器の能力範囲(流量 x 圧力)で表現できますが、 理解しやすいのはグラフ形式です(『23ページ図10』)。図からわかるよ うに、1290 Infinity ポンプの能力範囲は、1200 bar での動作と (2 ml/min 流量以下の場合)、800 bar への減圧(流量を 5 ml/min に増 加させた場合)です。この範囲は市場に出回っているすべての既存の UHPLC システムの動作範囲をカバーしており、これらのシステムから 1290 Infinity システムへメソッドを直接移管することを可能にしています。



2 Agilent 1290 Infinity LC システムの製品説明

Agilent 1290 Infinity LC システムの新機能

圧力の範囲を拡大したことによって、最新のサブ 2 マイクロンの粒子がサ ポートできるようになりました。そのため、長いカラムでは高い分解能、 短いカラムでは流量の増加に伴う高速分離を獲得できます。流量の範囲の 拡大によって、流量が多い場合に、従来のメソッドだけでなく、表面が多 孔質(または薄膜)の充填剤(たとえば、Poroshell)でも使用できるよ うになりました。この種の充填剤は最近では、効率的な分離用の STM 剤へ の代替アプローチとして着目されるようになってきています。流量の範囲 が拡大したことによって、分離に最適の内径を持つカラムを選択できるよ うになりました。すなわち、特定の MS システムの要件である低流量アプ リケーション用に 2 mm の内径か、以前からある LC 用またはローダビリ ティを増加させるために最大で 5 mm の内径 (通常は、4.6 mm)を選択で きます。流量の範囲が拡大したことで、最新の研究成果に対応できるよう になりました。その際には、グラジエント分離の効率を向上させるために 最大流量が使用できるという利点が活かされます(『Petersson et al., J. Sep. Sci, 31, 2346-2357, 2008, Maximizing peak capacity and separation speed in liquid chromatography』を参照)。

新しいダイオード アレイ検出器は、革新的なセル設計による優れたベース ライン特性と使い勝手と同時に、従来にはなかった優れた感度を備えてい ます。

オートサンプラでは、高圧力、低ボリュームアプリケーション用にアップ グレードされた、低キャリーオーバの可変ボリューム注入のために、定評 あるアジレントのフロースルー設計が採用されています。完全に新しいモ ジュールであるフレキシブルキューブは、オートサンプラに追加すること によって、固定ループ注入機能に対して、ディレイボリュームを最小化し、 ニードルシートのバック フラッシュのような性能改善を行います。

システムコンポーネント

Agilent 1290 Infinity バイナリポンプ

Agilent 1290 Infinity バイナリポンプ バルブクラスタ コントロールで は、LC 溶媒を非常に高い圧力と多くの流量で送り込む際の問題点を解決す るために、次のような新しいテクノロジが採用されています。すなわち、 過酷な使用に耐えるピストン上のドライブのモータ、負荷に耐えるだけで なくシールの放熱を促進するピストン用の新素材、マイクロフルイディッ ク熱交換器、マイクロ流体ミキサー Jet Weaver などです。このポンプは、 0.05 - 5 mL/min の範囲の流量をサポートします(最大圧力は、 1200 bar)。

Agilent 1290 Infinity バイナリポンプ バルブクラスタ コントロールモ ジュールには、2 つの同等な高圧 (1200 bar) ポンプ、2 チャンネルのデ ガッサと 4 チャンネルの注入溶媒切り替えバルブ、単一ハウジングに統合 された自動パージバルブと低ボリュームミキサー Jet Weaver が含まれま す。デガッサによって、流量が安定 (特に低流量時)し、検出器の感度が 向上します。



図 11 Agilent 1290 Infinity バイナリポンプ

Agilent 1290 Infinity LC システムマニュアルおよびクイックリファレンス

2 Agilent 1290 Infinity LC システムの製品説明

システムコンポーネント



図 12 1290 Infinity バイナリポンプの部品名と概略図

各ポンプヘッドは、新しいファームウェアコントロールと、ポンプの放熱 効果がある新しいピストン素材であるシリコンカーバイドを使った、デュ アルプランジャ直列型設計です。プライマリピストンをセカンダリピスト ンに結合しているキャピラリは、高い圧力と多くの流量で生成される熱を 放熱するための統合熱交換器を持っています。各ポンプヘッドは、プライ マリピストンチャンバに、パッシブインレットバルブ 1 つと、パッシブア ウトレットバルブ 1 つを持っています。各ピストンは、65000 ステップ のモーターで個別かつ正確に駆動され、各ステップで 300 ピコリッター の入れ換えを行います。

ピストンの移動は、圧力の脈動のアクティブダンピングがリップルフリー フローになるように、フィードバックループでインテリジェントに制御さ れます。ピストンドライブは、脈流のない状態を保つために、溶媒の圧縮 率特性とシステムの水力学特性を自動的に調整します。このような機能や、 ピストンの移動で発生する圧力の脈動を緩和するための円滑動作コント ロール機能、そして少ない容量の効率的なミキシングによって、UV シグナ ルのポンプノイズが最小限になることが保証されます。ポンプ内の専用マ イクロプロセッサが、静的/動的パラメータに基づいて、円滑な動作コン

Agilent 1290 Infinity LC システムの製品説明 2

システムコンポーネント

トロールとピストン動作の最適化をリアルタイムで制御します。クロマト グラフの性能の他に、これらの機能があることで、ポンプは非常に静かに 動作します。

移動相で高濃度の緩衝液を使用する場合は、ポンプシールの寿命を延ばす ためのアクティブシールウォッシュオプションを利用できます。

溶媒切り替えバルブを使えば、各チャンネルで 2 つの溶媒のうちの 1 つ を選択して、バイナリ混合物(アイソクラティックまたはグラジエント) を組成できます。バイナリグラジエントは、パージバルブでポンプ A とポ ンプ B の溶媒を高圧で混合することで生成されます。ソフトウェア制御の もとでパージバルブを使えば、ポンプヘッドから新しい溶媒をパージする ために、フローを廃液系に切り換えられます。システムの圧力を監視する ために、パージバルブにはセンサーが付属しています。



図 13 Jet Weaver ミキサー

ポンプの流路は、グラジエントディレイを最小化するように最適化されて おり、マイクロ流体テクノロジを用いた革新的なミキシングシステムが採 用されています。Jet Weaver として知られるミキサーでは、フローを完全 に混合できるようにするため、多層マイクロマイクロフルイディックチャ ンネル (120 μm x 120 μm)のネットワークが使われています。Jet Weaver には、2 つの標準ボリュームがあります。35 μ1: 通常の UV 検出アプリ ケーション用。100 μ1: UV 検出で TFA を使うときのように条件が厳しい 場合。これらの要求度の高いアプリケーションでベースラインのノイズを 徹底的に低減するために、380 μL の Jet Weaver も用意されています。MS 検出の場合は、多くの場合に Jet Weaver は不要です。ポンプ流路のベー ス 10 μl ボリュームしか使わなくても、十分に混合できます。このポンプ

2 Agilent 1290 Infinity LC システムの製品説明 システムコンポーネント

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。

1290 Infinity バイナリポンプには追加バルブレールが付属しており、ポ ンプの右側または左側に取り付けることができます。これらのバルブレー ルには、追加 12 段階溶媒切り替えバルブを最大で 2 つ取り付けられま す。そのため、バイナリグラジエントの分析メソッド開発では、最大で 26 種類の溶媒を使うことができます。外部溶媒切り替えバルブをポンプの ユーザーインタフェースに組み入れて、溶媒を簡単に名前で選択できるよ うにするための特殊な「クラスタドライバ」が用意されています。

28

1290 Infinity クォータナリポンプ

一方で、1290 Infinity クォータナリポンプは、ポンプヘッドを 1 つだけ 装備し、プログラムされたグラジエントに従って溶離液を分割するための 追加のマルチチャンネル グラジエントバルブ (MCGV)を備えています。こ の低圧混合の原理に従って溶媒がインレットウィーバーに集まるため、溶 媒はポンプヘッドの前およびポンプヘッド内で既に混合されています。

ポンプヘッドは、1290 Infininty バイナリ ポンプのものと同じであるため、同じパフォーマンス値を有し、技術的にも細部にわたり同等です。また、アクティブシールウォッシュも取り付けられているため、高濃度の緩 衝液の使用時でもポンプシールの寿命が延長されます。

分析中には、4 チャンネルのストリームセレクタである多目的バルブに溶 媒が入る前に圧力センサが圧力をモニタします。このバルブは、下図に示 すさまざまな機能を可能にしています。多目的バルブは、分析中に常時使 用されるインラインフィルタを備え、また取り付けが簡単で溶媒の最適な 混合を確保するオプションの 380 µL Jet Weaver ミキサー、オプションの リストリクションキャピラリ、および排液システム接続部を取り付けるこ とができます。

2 Agilent 1290 Infinity LC システムの製品説明 システムコンポーネント

表 2 バルブの機能 ポンプ サンプラ す う て イルタ



図 14 標準アプリケーション

図 15 エクストラミキシングボリューム のセットアップ





図 17 インラインフィルタのバック フ ラッシュ

図 16 自動パージ機能

Agilent 1290 Infinity LC システムの製品説明 2

システムコンポーネント

標準アプリケーション(1)はほとんどの分析に使用されます。一方で、エ クストラミキシングボリュームのセットアップ(2)は、Agilent Jet Weaver ミキサーの使用によって混合パフォーマンスと依存性 UV ベースラ インを大幅に強化できるような、ベースラインが大きな意味を持つアプリ ケーションに適用されます。自動パージ機能(3)とインラインフィルタの バック フラッシュ(4)の構成も可能です。これによりフィルタを清潔に 保つことができ、寿命が延長されます。

2 Agilent 1290 Infinity LC システムの製品説明 システムコンポーネント

Agilent 1290 Infinity オートサンプラ

Agilent 1290 Infinity オートサンプラでは、可変ボリューム注入用に定 評あるアジレントのフロースルー設計が採用され、画期的な性能が実現で きています。メタリングデバイスのシールとニードルシート内で新しい不 活性剤を使うと、極めて低いキャリーオーバを達成できます。流路の水圧 低下機能はグラジエントを高速化することに貢献し、オーバーラップ注入 機能と自動バルブ切り替え (ADVR)機能は、サイクルタイムの短縮と、グ ラジエントのカラムへの高速送液に貢献します。このシステムは、サンプ ル溶液のセットボリュームを無駄なく正確に吸引し、サブマイクロリッ ターから最大注入量 40 µl の範囲にわたって高い再現性を実現します。標 準で取り付けられている注入キャピラリを使えば、20 µl だけ注入できま す。



図 18 1290 Infinity オートサンプラの注入ステップ概略図

Agilent 1290 Infinity LC システムマニュアルおよびクイックリファレンス

Agilent 1290 Infinity LC システムの製品説明 2

システムコンポーネント

完全に新規のオプションアドオンモジュールであるフレキシブルキューブ は、オートサンプラとシームレスに協調して、追加機能を提供します。新 規のフレキシブルキューブモジュールを追加する(500 µl シリンジポンプ を構成する)ことによって、1 つの低圧バルブと 2 つの高圧切り替えバル ブを構成できます。たとえば、シリンジポンプとフレキシブルキューブの バルブ操作を使ってサンプルループを構成すれば、フロースルー注入シス テムを固定注入ループで置き換えられます。この構成の利点は、オートサ ンプラからディレイボリュームを除去できることです。つまり、高速グラ ジエント、ハイ スループットの状況に適した構成となります。トレードオ フは、可変ボリューム注入の自由度が消滅し、サンプルによってはループ 内のフラッシングでは、注入ニードルシートは、注入後にフレキシブル キューブを使って、汚れたサンプルが原因となる複雑な化合物のキャリー オーバや詰まりを防止します。

オートサンプラのサンプルラックには、2 ml バイアルと 2 つの取り外し 可能なトレイ用の 10 ポジションがあります。2 つの取り外し可能なトレ イは、それぞれ独立に以下の用途で使用できます。

- 54 ポジションを持つ 2 ml バイアル
- 96 ウェルポジションを持つマイクロプレート(高さは可変)
- 384 ウェルポジションを持つマイクロプレート(高さは可変)

必要に応じてオートサンプラにオートサンプラ温度コントロールモジュー ルを追加すれば、4 °C~40 °Cの範囲で温度調節ができるようになりま す。

Agilent 1290 Infinity カラム コンパートメント

Agilent 1290 Infinity カラム コンパートメント (TCC) は、周囲温度が -10°C から、100°C (2.5 ml/min の場合)または 80°C (5 ml/min の 場合)までの範囲のときに、温度が制御できます。温度安定性の仕様は ± 0.05°C であり、精度仕様は ± 0.5°C (キャリブレーション時)で $_{\sigma}^{1}$.この条件は、サーモスタットベインの接触部の伝導、カラム環境内 の静止空気温度、そして最も重要ですが、移動相を、カラムに挿入する前 に熱交換器内をくぐらせることによって、前もって加熱(または冷却)し ておくことを組み合わせれば達成できます。各 TCC 内には、300 mm まで の長いカラム用の温度帯と、100 mm 以下の短いカラム用の温度帯の 2 つ の独立した温度帯があります。

このモジュールには、1.6 μ l の低拡散熱交換器が付属しています。各バル ブキットは、各カラム用の追加低拡散熱交換器を持っています。低拡散熱 交換器は、TCC 内に最大 4 台まで自由に取り付けることができます。従来 の HPLC 操作を行う場合は、3 μ l と 6 μ l の内蔵熱交換器を使うこともで きます。

各 TCC には、1 つの内部バルブドライブを収容できます。バルブ切り替え アプリケーションはこのドライブを使って、2 つのカラムを単純に切り替 える自動的なカラム再生、サンプルの準備、カラムバックフラッシングな どを実行します。各バルブヘッドは、すべての必要なキャピラリ、追加低 拡散熱交換器、その他の部品を含む完全なキットとして提供されます。

切り替えバルブは、バルブに接続すれば、非常に使いやすく、柔軟性がある操作性を提供します。押し込めば、クイックチェインジバルブのドライブユニットが前方向にスライドし、簡単にアクセスできるようになります(『35ページ図19』の左図を参照)。異なるアプリケーションの場合は、ユーザーがドライブメカニズムの代替バルブヘッドを交換できます(『35ページ図19』の右図を参照)。バルブヘッドの上部にある RFID タグに注意してください。

¹ いずれの仕様も、周囲温度 25 ℃、設定温度 40 ℃、流量範囲 0.2 ~ 5 ml/分の蒸留 水に対して有効です。

Agilent 1290 Infinity LC システムの製品説明 2

システムコンポーネント



図 19 TCC 内のクイックチェインジバルブ

TCC は最大 3 台まで「クラスタ化」できます。この構成では、自動メソッ ド開発用に 8 つのカラムを切り換えるような高度なアプリケーションに対 応したり、異なるアプリケーションに追加カラムを提供することができま す。したがって、使用するカラムは、単純なメソッド パラメータになりま す。これには 2 つの 8 ポジション /9 ポートバルブヘッドが必要ですが、 これは 2 つの TCC につき、1 つしか付属していません。クラスタ構成の TCC は、操作が楽になるように、ソフトウェアによって、1 つのインタ フェースしか持たない 1 つのユニットとして表現されます。

以前の設計に比べて改善された点は、断熱効果が向上したこと、キャピラ リガイドが改善されたこと、そして「ドアオープン」センサーが追加され たことです。このセンサーによって、メソッドでドアが閉じられた状態を 定義できるようになりましたが、これは特に低温度 / 高温度アプリケー ションで役に立ちます。

2 Agilent 1290 Infinity LC システムの製品説明 システムコンポーネント

1290 Infinity ダイオードアレイ検出器

1290 Infinity ダイオードアレイ検出器は、光学流体導波路テクノロジを 採用したカートリッジセルを用いた新しい光学設計です。このテクノロジ は、標準または超高速 LC アプリケーションに対して、高感度、低拡散、 広い線形範囲、そして非常に安定したベースラインなどの属性を提供しま す。Agilent Max-Light カートリッジセルでは、コーティングされていな いフューズドシリカキャピラリに沿った全反射を利用することによって、 光の透過率が著しく改善されました。つまり、セルボリュームの拡散効果 を通じて、分離性能を犠牲にすることなく、従来よりもはるかに優れた感 度を達成しています。この設計により、屈折率や熱効果が原因となるベー スライン変動を最小化できるため、ピーク面積の積分の信頼性が向上しま す。



1024 素子ダイオードアレイ

図 20 1290 Infinity ダイオードアレイ検出器を経由する光パス

このモジュールには、周囲温度の影響への耐性を強化するための、電子温 度コントロールも採用されています。Max-Light カートリッジセルの水力 ボリュームは非常に小さいものの、光路長は標準の 10 mm です。ただし、 さらに高感度が必要な場合には、光路長が 60 mm の別の Agilent Max-Light 高感度セルも用意されています。セルは、セルホルダーの前後 にスライドさせることで簡単に交換でき、光学作業台内で自動調整されま す。DAD の光源は、重水素ランプです。動作波長範囲は、190 ~ 640 nm を カバーしています。これは、1024 ダイオードで構成されるダイオードアレ
Agilent 1290 Infinity LC システムの製品説明 2

システムコンポーネント

イで検出されます。分光器への入力は、1~8 nm のスペクトル分解能を持 っプログラム可能な光学スリットを経由して行われます。これは通常、範 囲の中位点で動作しますが、高スペクトル分解能のためには 1 nm まで狭 くすることができます。ただし、液相 UV スペクトルでこれが必要になる ことはめったにありません。また最大光透過率と最小シグナルノイズが必 要な場合には、8 nm にすることができます。

クロマトグラフ シグナルは、モジュール内のファームウェアによって、ダ イオードアレイのデータから抽出されます。最大で 8 つの独立シグナルを 定義できます。各シグナルは、シグナル波長、ダイオードバンチングバン ド幅、そして、必要に応じて、リファレンス波長とバンド幅で定義します。 シグナルの出力で、最高速(最狭)クロマトグラフ ピークを正確に記録 できるのは、最高で 160 Hz (160 データ数/秒)の場合です。また、この モジュールは、同じレート 160 Hz で全範囲のスペクトルをデータシステ ムに出力できます。

規制対象下のラボでは、すべてのメソッド パラメータを記録しておく必要 があります。1290 Infinity DAD は、機器の設定値だけでなく、ランプや フローセルカートリッジ内の RFID タグ (無線周波数識別タグ) も記録し ます。そのため、システムには、これらの重要なコンポーネントの ID や 変数も記録されます。

2 Agilent 1290 Infinity LC システムの製品説明 システムコンポーネント



3

Agilent 1290 Infinity LC システムマニュアルおよびクイック リファレンス

Agilent 1290 Infinity LC システ ムの最適化

ディレイボリュームとエクストラカラムボリューム 40 ディレイボリューム 40 エクストラカラムボリューム 41 最適ディレイボリュームの設定方法 42 注入量を増加させる方法 51 ハイスループットを達成する方法 53 高分解能を達成する方法 56 感度を向上させる方法 59 キャリーオーバを最少にする方法 67 カラムの詰まりを回避する方法 69

この章では、理論の適用方法と、LC システムの特徴を活かして最適分離を達成する方法について説明します。



3 Agilent 1290 Infinity LC システムの最適化 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、カラムへの到達時間が変化します。 ディレイは、流量とシステムのディレイボリュームに依存します。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設定を使ってレ 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、低流量と小さ 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の転送 に対しては、最大の影響を与えます。したがって、高速グラジエント分離 の場合は、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 分析検出でしばしば使われる内径が狭いカラム(たとえば、2.1 mm の内径)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。

たとえば、5 µm の充填剤を使う HPLC メソッドでは、流量 1 ml/分 が、 通常、内径 4.6 mm のカラムで使われ、流量 0.2 ml/分 が、内径 2.1 mm のカラムで使われます (カラム内の線速度が同じ場合)。一般的な 1000 µ1 のディレイボリュームを持ち、2.1 mm のカラムを使うシステムに は、最初から 5 分 の「隠れた」アイソクラティックセグメントがありま す。一方、600 µ1 のディレイボリュームでは、ディレイは 3 分になりま す。このようなディレイボリュームは、1 分または 2 分程度の分析時間に 対しては大きすぎます。サブ 2 マイクロン充填剤を使う場合は、最適の流 量 (Van Deemter 曲線による)が若干増加し、高速クロマトグラフィでは この流量の 3~5 倍の流量を使用できるので、ディレイ時間は 1 分ほど

ディレイボリュームとエクストラカラムボリューム

で済むようになります。ただし、ディレイ時間を、予定している分析時間 の数分の1にするには、ディレイボリュームをさらに減らす必要がありま す。Agilent 1290 Infinity LC システムでは、ポンプの流路のディレイボ リュームが少なくなったこと、Jet Weaver ミキサーのボリュームが少いこ と、オートサンプラを経由する流路のボリュームが少なくなったことで、 これが実現できています。

エクストラカラムボリューム

エクストラカラムボリュームは、分離分解能を低下させ、それ故に、最小 化する必要があるピーク拡散の原因です。ピーク拡散を最小限に抑えるに は、カラムの内径を小さくすることに比例して、エクストラカラムボ リュームも小さくする必要があります。

液体クロマトグラフでは、エクストラカラムボリュームは、オートサンプ ラ、カラム、検出器間の接続チューブに依存します。検出器内では、フ ローセルのボリュームに依存します。Agilent 1290 Infinity LC システム では、チューブ接続の内径を狭くした(0.12 mm)こと、カラムコンパー トメント内の熱交換器と検出器内の Max-Light カートリッジセルのボ リュームを小さくしたことによって、エクストラカラムボリュームが最小 となっています。

3 Agilent 1290 Infinity LC システムの最適化 長海ディレイボリュームの設定支法

最適ディレイボリュームの設定方法

最適ディレイボリュームの設定方法

『42ページ 表 3』と『43ページ 表 4』に、Agilent 1290 Infinity LC シ ステム内のシステムディレイボリュームの原因となるコンポーネントボ リュームを示します。Jet Weaver ミキサーが付属する Agilent 1290 Infinity バイナリポンプ バルブクラスタ コントロール、1290 Infinity オートサンプラ、カラムコンパートメントを使用する標準構成では、シス テムのディレイボリュームは、およそです 125 µl です。この標準的な ディレイボリュームは、大多数のアプリケーションには適しています。た とえば、サブ 2 マイクロン粒子が充填された 50 mm x 2.1 mm のカラムと 中間的な流量 0.6 ml/min で高速分離を行った場合は、約 0.2 min の一般 的なグラジエント時間が生成されますが、この値は、2~3 分のグラジエ ント時間の場合は一般的には許容範囲です(『44ページ 表 6』を参照)。 流量をカラムボリュームを用いて考えることが役立つ場合があります。た とえば、『44ページ 表 7』の 0.6 ml/min カラムでは、システムで 1 分当 たり 6 カラムボリュームが流れており、ディレイボリュームはカラムボ リュームの約 1.2 倍であることが観察できます。

Agilent 1290 Infinity クォータナリポンプ、1290 Infinity オートサン プラ、およびカラム コンパートメントを使用する構成のディレイボリュー ムは 430 µL であり、結果としてディレイタイムは 0.7 min となります。 この時間は、グラジエント時間が 3 min の場合の許容限界値です。

コンポーネント	ディレイボリューム (μL)
バイナリポンプ	10
Jet Weaver ミキサー (標準)	35
バイナリポンプ + Jet Weaver	45
クォータナリポンプ	350
クォータナリポンプ + V380 Jet Weaver	500
オートサンプラ(固定 5 µl ループ)	5
オートサンプラ(標準、可変ボリューム)	80

表 3 1290 Infinity LC モジュールのディレイボリューム

Agilent 1290 Infinity LC システムマニュアルおよびクイックリファレンス

最適ディレイボリュームの設定方法

コンポーネント	ディレイボリューム (μL)
カラムコンパートメントの低拡散熱交換器	1.6
接続チューブ(0.12 mm 内径)100 mm につき	1.1

表 3 1290 Infinity LC モジュールのディレイボリューム

表 4 1290 Infinity バイナリ LC システム構成のディレイボリューム

システム構成 ¹	ディレイボリューム (μL)
バイナリポンプ + 固定ループオートサンプラ(MS のみ)	20
バイナリポンプ + Jet Weaver + 固定ループ	55
バイナリポンプ + 標準オートサンプラ(MS のみ)	90
バイナリポンプ + Jet Weaver + オートサンプラ	125

¹ システム構成内の接続のために 5 µl を追加

表 5 1290 Infinity クォータナリ LC システム構成のディレイボ リューム

システム構成 ¹	ディレイボリューム (μL)
クォータナリポンプ + 固定ループオートサンプラ (MS のみ)	360
クォータナリポンプ + 標準オートサンプラ (MS の み)	430
クォータナリポンプ + V380 Jet Weaver + 固定 ループ	510
クォータナリポンプ + V380 Jet Weaver + オート サンプラ	580
1	

¹ システム構成内の接続のために 5 µl を追加

最適ディレイボリュームの設定方法

流量 (m1/分)		システム	ディレイ	'ボリュ-	-4 (マ	イクロリ	ッター)	
	20	55	90	125	360	395	430	465
			ア	イレイタ	74(3)	子)		
0.2	0.10	0.28	0.43	0.60	1.80	2.15	2.55	2.90
0.4	0.05	0.14	0.21	0.30	0.90	1.08	1.28	1.45
0.6	0.03	0.09	0.14	0.20	0.60	0.72	0.85	0.97
0.8	0.03	0.07	0.11	0.15	0.45	0.54	0.64	0.73
1.0	0.02	0.06	0.09	0.12	0.36	0.43	0.51	0.58
1.5	0.01	0.04	0.06	0.08	0.24	0.29	0.34	0.39
2.0	0.01	0.03	0.04	0.06	0.18	0.22	0.26	0.29
3.0	0.01	0.02	0.03	0.04	0.12	0.14	0.17	0.19
4.0	0.01	0.01	0.02	0.03	0.09	0.11	0.13	0.15
5.0	0.00	0.01	0.02	0.02	0.07	0.09	0.10	0.12

表 6 グラジエントがカラムヘッドに到達するまでのシステムディレイ タイム

表 7 典型的な寸法のカラム内の概算液体ボリューム (多孔率 = 0.6 を仮定)

カラム内径 (mm)	カラムの長さ (mm)				
	30	50	100	150	250
	カラゴ	ムボリューム	- 液体相(-	マイクロリッ	ター)
2.1	62	104	208	312	520
3.0	127	212	424	636	1060

Agilent 1290 Infinity LC システムマニュアルおよびクイックリファレンス

最適ディレイボリュームの設定方法

カラム内径 (mm)		力	ラムの長さ(mm)	
4.0	226	377	754	1131	1885
4.6	299	499	997	1496	2493

表 7 典型的な寸法のカラム内の概算液体ボリューム (多孔率 = 0.6 を仮定)

Agilent 1290 Infinity バイナリ LC システムのみが対応可能な 0.5 min 程度の非常に早いグラジエント分析を実施するために、システムの物理構 成を変更することなく、システムのディレイボリュームを容易に縮小でき ます。縮小は、オートサンプラの動作を変更するだけで済みます。

Agilent 1290 Infinity オートサンプラの 80 ul ディレイボリュームは、 メタリングデバイスを経由する注入バルブからの流路、ニードル、ニード ルシート、注入バルブに戻る接続キャピラリに依存します(『32ページ 図 18』を参照)。注入を行うためにバルブはメインパスからバイパスへ切 り替えるため、メタリングデバイスはニードルキャピラリにサンプルを吸 引できるようになります。バルブがメインパスに切り替わるときに注入が 行われ、サンプルがカラム内にフラッシュされます。解析中にバルブはこ のポジションに留まり、オートサンプラはフラッシュされ続るため、グラ ジエントがカラムに到達するにはこのディレイボリューム内を流れなけれ ばなりません。これは、注入が完了して、注入されたサンプルがカラムに フラッシュされた後で注入バルブをメインパスからバイパスへ切り替える ことで回避できます。実際に切り替わるのは、注入が終わって数秒経過し た後です。この機能は、オートサンプラの設定メニューで、[自動バルブ 切り替え] (ADVR) 機能を選択すればアクティブになります。フラッシュア ウト係数(通常は、注入量の5倍)は、バイパスへ切り替える前にイン ジェクタからサンプルをフラッシュするための十分な時間を確保するため のパラメータです。この係数によって、システムディレイボリュームは、 125 µl から 50 µl に減少します。

ADVR を使う場合は、注入が行われる時点で、ポンプでグラジエントが開始 されている必要があります。グラジエントがオートサンプラに到達済みか どうかも問題になります。到達済みの場合は、グラジエント内に小さなス テップが生じるためです。これはディレイボリュームがフラッシュアウト ボリュームよりも少ない場合に発生します。これが常に問題となるわけで はありませんが、メソッド移管時には考慮が必要です。フラッシュアウト 係数が 5 で、注入量が 10 µl の場合は、オートサンプラはバイパスに切

Agilent 1290 Infinity LC システムマニュアルおよびクイックリファレンス

45

最適ディレイボリュームの設定方法

り替える前に、50 µ1 を通過させることになります。ディレイボリューム 50 µ1 は、グラジエントが注入バルブに到達したばかりであることを意味 します。注入量が少なければ影響はありませんが、多ければグラジエント 内に小さなステップが生じることになります。使用中の流量は、ADVR の使 用可否にも影響します。ディレイタイムが 0.2 ml/min の場合は、短縮さ れるディレイタイムは 21 秒ですが、1.0 ml/min の場合は 4 秒になりま す。

ADVR 機能は、キャリーオーバ問題を引き起こすことがわかっている化合物 を含むアプリケーションには適していません。

カラム コンパートメントでピーク拡散とディレイボリュームを最小化する には、低拡散熱交換器を取り付けておく必要があります。低拡散熱交換器 はキャピラリキットに含まれており、低拡散アプリケーションでの使用が 推奨されています。通常のキャピラリキットには、内径が 0.12 mm の狭い キャピラリも含まれています。内蔵の 3 µl と 6 µl の熱交換器は従来の 機器との互換性のためだけに用意されています。したがって、使用するの は、汎用メソッドを実行する必要がある場合だけにしてください。ただし、 汎用メソッドを実行する場合でも、低拡散熱交換器を使うことは可能です。

Agilent 1290 Infinity ダイオードアレイ検出器の分解能を劣化させない ために、Max-Light カートリッジセルには低拡散ボリューム (s ボリュー ム 1.0 μ l) が付属しています。そのため、ボリュームの最適化は不要で す。高感度が必要なために代替の Agilent Max-Light 高感度セルを使う場 合は、セルボリュームは 3 mm と 4.6 mm の内径のカラム用に最適化する 必要があります。

ポンプが正常に動作するためには、ポンプのセットアップ画面で正しい溶 媒を設定する必要があります。インテリジェントコントロールによって、 圧力リップルは自動的に最小になるように調整されますが、溶媒圧縮率は 高圧時の正しい流量を極めて正確に維持することに影響を与えます。これ により、使用中の移動相に対して常に正しい圧縮率値を適用することが保 証されます。キャリブレーション機能は、Agilent 1290 Infinity バイナ リポンプ バルブクラスタ コントロールおよびクォータナリポンプで使用 できます。

1290 Infinity バイナリ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 ムから 35 µl を削除することによって、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。ディ

最適ディレイボリュームの設定方法

レイタイムがシステムに与える影響については、『44 ページ 表 6』を参照 してください。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。Jet Weaver を 外す場合は、バッファやその他の添加剤が残らないように、溶媒でフラッ シュしてください。

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。Jet Weaver ミキサーには、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 µl から高ボリューム 100 µl への切り換えは、インストール解除、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。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 µl 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。UV ベースラインノイズを可能な限り低減させる必要のある要求度の高いアプリケーションには、標準の Jet Weaver ミキサーの代わりに取り付けられる 380 µL Jet Weaver ミキサーを使用できます。

1290 Infinity バイナリポンプの Jet Weaver の交換手順を、『「1290 Infinity バイナリポンプの Jet Weaver の交換」48 ページ 図』に示します。

Agilent 1290 Infinity クォータナリポンプは構成と混合原理が異なるため、物理的なディレイボリュームがはるかに大きくなり、標準アプリケーションで Jet Weaver ミキサーを追加する必要はありません。ただし、TFA アプリケーションなどのベースラインが重要な意味を持つアプリケーションの場合は、オプションの 380 µL Jet Weaver ミキサーを取り付けることも可能です。オプションの Jet Weaver ミキサーには、1290 Infinity クォータナリポンプの設計に適合させるための異なるハウジングが採用されています。

1290 Infinity クォータナリポンプの Jet Weaver の取り付け手順を、 『「1290 Infinity クォータナリポンプの V380 Jet Weaver の取り付け」49 ページ 図』に示します。

1290 Infinity バイナリポンプの Jet Weaver の交換



1290 Infinity クォータナリポンプの V380 Jet Weaver の 取り付け



Agilent 1290 Infinity LC システムマニュアルおよびクイックリファレンス

最適ディレイボリュームの設定方法



注入量を増加させる方法

注入量を増加させる方法

Agilent 1290 Infinity オートサンプラの標準構成には、最大で 20 µ1 の 注入用の可変ボリュームサンプルループが含まれています。メタリングデ バイスは、最大で 40 µ1 のボリュームを注入できます。これを利用するた めに、サンプルループカートリッジを交換できます(詳細は、1290 Infinity オートサンプラマニュアルを参照)。オートサンプラで発生する システムディレイボリュームは、それに応じて増加します。

メソッドを大きなカラムから小さなカラムにスケールダウンする場合は、 メソッドの性能を維持するために、メソッド変換でカラムのボリュームに 比例した注入量を減らす必要があります。これによって、注入量の割合を カラムと同じ割合に保ちます。この操作は、注入する溶媒が開始移動相よ り強い(有機溶媒が多い)場合は特に重要です。多いと、必ず分離性能に 影響があります(特に、溶出時間の短い分析ピークの分離性能)。場合に よっては、これがピーク拡散の原因になります。一般に、注入する溶媒を 最初のグラジエント混合と同じか、それよりも弱くしておく必要がありま す。これは、注入量を増加できるかどうか、また増加できる場合は、どの 程度増加できるのかといったことに関連があります。注入量の増加を試み る場合は、拡散が増加している徴候(ピークの広がりまたは歪みやピーク 解像度の減衰)がないかどうかを確認する必要があります。弱い溶媒に注 入すると、グラジエント開始時にカラムのヘッドにある対象化合物を濃縮 する結果となります。逆に、開始移動相よりも強い溶媒に注入すると、増 加した注入量が、グラジエントよりも前にあるカラムに沿って対象化合物 を分散させるため、ピーク拡散と分解能の低下を招きます。

したがって、注入量を決定する際に最も考慮する必要があることはカラムの内径です。これがピーク拡散に最も影響があるからです。狭いカラムでは、ピーク拡散が減少するため、広いカラムで多くの注入を使った場合よりもピークの高さが高くなります。内径が2.1 mmのカラムでは、一般的な注入量の範囲は 5~10 µl ですが、前述のように、これは対象化合物の化学的性質と移動相に大きく依存しています。最大注入量のおよその目安は、カラムのボリュームで判定できます(『』を参照)。最適の分離性能とピーク拡散を保つには、カラムボリュームの約5%のグラジエント分離注入量が必要になります。

試料量を増加させるための 1 つの方法は、切り替えバルブで選択されるト ラッピングカラムを使用して、注入物を分析カラムに切り替える(注入す

注入量を増加させる方法

る)前に、注入物を取得してそれを濃縮することです(『』を参照)。バ ルブは、温度調節機能付きカラムコンパートメントの便利な位置に取り付 けられます。



ハイスループットを達成する方法

ハイスループットを達成する方法

数百または数千の注入が必要となるハイスループット(HT)環境で運用しているラボがあります。このような状況では、サイクルタイムの短縮が最大の課題であり、1回の注入で、たとえ数秒しか短縮できないとしても、作業の完了までに必要な合計時間は、著しく、そして有意義に減少させます。高速サイクルタイムとハイスループットオペレーションを達成するための、重要なステップは次のとおりです。

- 高速分離を利用する
- 注入をオーバラップさせる
- 平衡時間を最小化する
- カラム再生を交互に行う

ハイスループットオペレーションを達成するための最初のステップは、サ イクルタイムの短い高速クロマトグラフィメソッドを使うことです。この 目的のためには、粒子径が 1.8 µm の充填剤を使う短いカラムを使うこと が理想的です。短いカラムは高効率だからです。メソッドがアイソクラ ティック分離を含んでいる場合は、新しい分析に移るたびに必要となる平 衡時間が生じないため、短いカラムを使うことによって、高速サイクルタ イムが実現できます。ただし、多くの場合、サンプルの範囲や複雑さのた めに、グラジエントメソッドが使われます。メソッドの開発中は、グラジ エント範囲は、分離を行うのに必要な最小限に保つ必要があります。多く の「オープンアクセス」システムでは、どのような不明な化合物にでも対 応することが必要なことから、5 % ~ 95 % の有機溶媒を用いるグラジエ ントが使われます。ハイスループット環境では、予想される化合物の範囲 を狭めることができるか検討する必要があります。これは分析時間を短縮 するだけでなく、新しい分析に移るたびに必要となる平衡時間も短縮する ことに貢献するからです。

サイクルタイムは、次のようにいくつかの要素で構成されます。サイクル タイム = 注入 + 分離 + 平衡化 + データ処理

処理対象のサンプル数が大きい場合は、サイクルタイムを僅かに短縮した だけでも、作業完了までの合計時間の大きな低減につながります。そのた め、データ処理はオフラインで行って、システムはサンプルの分析とデー タ収集に専念できるようにします。

ハイスループットを達成する方法

サンプルの取り出しを急ぐと再現性を損なうことさえ忘れなければ、注入 の速度は最適化できます。多くの場合に、使用するサンプル量が、範囲の 下限となっている傾向があるため、限界ゲインが求められます。注入時間 の大部分は、ニードルをバイアルとフラッシュポート間で移動させるのに 必要な時間です。このような操作は、先行する分離を行っている最中に実 行できます。これは「オーバラップ注入」として知られており、 ChemStation コントロール ソフトウェアのオートサンプラ設定画面で有効 にできます。オートサンプラには、注入が終わったら、バイパスするため にフローの切り替えを指示できます。そして、たとえば、4 分間の分析で は、3 分後に次のサンプルを吸引して注入を準備するプロセスを開始でき ます。これにより、一回の注入で、通常は、0.5~1 分が短縮できます。 粘着性のある化合物の場合は、カラム平衡化の最中に行うことをお勧めし ます。これは、オートサンプラが次のグラジエント分析の開始条件を認識 するときと一致します。

カラム平衡化ステップは、サイクルタイムの大きな割合を占めます。通常、 次回の注入を準備するためにカラムボリュームを安定化させるまで、カラ ムでは 3~5 回のフラッシングが必要です。この操作はアプリケーション によっては、分離時間の 50 % 以上を占めます。これは本質的なプロセス ですが、自動交互カラム再生機能を使うことでサイクルタイムから除外で きます。そのためには、カラムコンパートメント内に 2 ポジション 10 ポート、1200 bar のバルブヘッド、それと同等の 2 番目の分析カラムが必 要です。また、2番目のポンプも必要です。1つのカラムが分離分析で使 われているときに、他のカラムは移動相グラジエントの初期混合でフラッ シュされます。そして、次の注入を開始するために、新たに平衡化された カラムが分析流路内に組み込まれます。2 つのカラムはこのように交互に 使われて注入シーケンスを構成します。2番目のポンプが必要になるのは、 カラム内のアイソクラティック混合物をフラッシュする場合だけです。そ のため、1290 Infinity ポンプよりも簡単な機構のポンプで済みます。た とえば、1200 Series アイソクラティックポンプで十分です。設定を、『55 ページ 図 22』に示します。

ハイスループットを達成する方法



Agilent 1290 Infinity LC システムマニュアルおよびクイックリファレンス

高分解能を達成する方法

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。この節では、以下の要素を検討することによって、分解能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、高速フロー
- 最小限のエクストラカラムボリューム
- ・ 注入溶媒とボリュームの最適化
- 十分に高速なデータ収集
- 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 = 分解能
- N = 段数(カラム効率の目安)、
- a = 選択性(2 つのピーク間)
- k₂=2 番目のピークの保留係数(旧称、キャパシティ係数)

分解能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(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で 評価できます。この節では、任意の固定相および移動相の分解能を向上さ せる方法について説明します。相の判断で自動メソッド開発システムを使

高分解能を達成する方法

うと、スカウティングの各ステップ内の分析を高速化するために、多くの 場合は短いカラムが選択されます。

分解能方程式によれば、次に重要な項は段数、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 1200 bar までの圧力とサブ 2 マイクロンの粒子に対応し、カ ラムの長さを 100 mm または 150 mm に延ばせるように設計された理由で す。また、100 mm と 150 mm のカラムを継ぎ足して、250 mm の長さのカ ラムを構成することもできます。分解能は N の平方根で増加するので、カ ラムの長さを倍にすれば、分解能は 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か らです。メタノール混合物は、アセトニトリル混合物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、良く使われますが、選択性はメタノールの方が優れて います(特に、500 Da 以下の小さな分子の場合)。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を変化させる可能性があることに 注意してください。経験すれば、選択性の増加または減少のいずれの原因 になるかがわかります。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、拡散を若干増加させ、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、分解能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 カラム内の最適の流量は、大きな粒子に対して は多く、流量が増加するに連れて平準化していくことを示しています。STM カラムの最適流量は、通常、以下のとおりです。2 ml/分(内径が 4.6 mm の場合)および 0.4 ml/分(内径が 2.1 mm の場合)。

アイソクラティック分離では、保留係数 k を増加させると、溶液の保留時 間が増加するため分解能が向上します。グラジエント分離では、保留は次 の方程式の 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

高分解能を達成する方法

- ・ $t_G =$ 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、ひいては分解能は、グラジエントを浅く(目安 としては、2~5%/分の変化)、流量を多く、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 \mathbb{C} 度 以上の温度で、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 を参照してください。)

エクストラカラムボリュームは、短くすれば拡散が減るので、分解能は向 上します。1290 Infinity LC システムの狭い内径(0.12 mm)のキャピラ リ(カラムと検出器間で最短の長さが使われていることを確認してくださ い)と Max-light カートリッジフローセルでは最適化済みです。

最後に、狭いピークを正確にプロファイルするのに十分な速度を持つデー タ収集であれば、分解能の向上に利することがあることを付記しておきま す。

58

感度を向上させる方法

分離メソッドの感度は、固定相 / 移動相の選択に依存しています。狭い ピークと安定したベースラインを持ち、ノイズが最小化された良好な分離 が目標となります。機器構成の選択も影響しますが、影響が最も大きいの は検出器の設定です。この節では、感度が以下の要素から受ける影響につ いて説明します。

- ・ ポンプミキサーボリューム
- 狭いカラム
- 検出器フローセル
- 検出器のパラメータ

また、検出器パラメータの説明でも、選択性と線形性についての関連ト ピックを説明します。

ポンプミキサーボリューム

UV 検出でベースラインノイズを最小限にするには、1290 Infinity ポンプ モジュールでは 35 µl の Jet Weaver を使う標準のディレイボリューム構 成をお勧めします。これは大部分のアプリケーションに適用できますが、 移動相で TFA を使用したり、多くの混合を必要とするような状況では、ミ キシングノイズを最小化するために、Jet Weaver ミキサーの大きなボ リュームサイドを使う必要があります。

カラム

感度は、シグナル対ノイズ (S/N) 比で表現されます。これは、ピークの高 さを最大化し、ベースラインノイズを最小化するために必要となるパラ メータです。ピーク拡散を減らせばピークの高さを維持できるので、エク ストラカラムボリュームは、短く狭い内径のキャピラリを使って、正しく 取り付けることにより、最小化する必要があります。小さな内径のカラム の使用は、ピークの高さを高くするので、限られたサンプル量しか使えな いアプリケーションには理想的です。同じサンプル量でも、小さな内径の カラムに注入した方が、カラムの内径が原因となる希釈が少なく、感度は 向上します。たとえば、カラムの内径を 4.6 mm から 2.1 mm に縮小すれ ば、カラム内の希釈がすくなくなるため、ピークの高さは理論的には

Agilent 1290 Infinity LC システムマニュアルおよびクイックリファレンス

59

感度を向上させる方法

4.7 倍 増加します。質量分析検出器の場合は、狭いカラムで流量を減らせば、イオン化の効率が上がるため感度が向上します。

検出器感度を向上させる方法

検出器には、性能を最適化するために使われるパラメータが多数あります。 以下のセクションでは、検出器のパラメータがパフォーマンス特性に与え る影響について説明します。

- フローセルが感度に与える影響
- 波長と帯域幅が感度、選択性、直線性に与える影響
- スリット幅が感度、スペクトル分解能、直線性に与える影響
- ピーク幅が感度と分解能に与える影響

フローセル

Max-Light カートリッジフローセルは、標準の 10 mm 光路長であり、最小ボリュームおよび分散用に最適化されています (s 容量 1.0 μ L)。これは、ノイズを最小にするための多光量転送機能を持っており、光学流体導波路が原因となるノイズを減らしています。これは、短く狭い内径のカラムから長く標準の内径(4.6 mm)のカラムまで、広い範囲の分析カラムで使用するのに適しています。一般的にピーク分散ボリューム(ピーク幅 x 流量で計算)は、このセルのボリュームに比べて、約 2 μ L 大きくなります(例えば 0.02 min x 200 μ L/min = 4 μ L)。

Max-Light 高感度セルは、60 mm の光路長を持っているため、アプリケー ションの条件に応じて、シグナル / ノイズの値は 3 ~ 5 倍増加します。 分散ボリュームは、標準セルに比べてわずかに増加します。

波長と帯域幅

検出器は、ダイオードアレイ検出を使って、190 nm ~ 640 nm の波長で吸 光度を一度に測定します。UV ランプは、全波長範囲にわたって高感度で す。ダイオードアレイ検出器 (DAD) は、検出時間ごとに最大で 8 つのク ロマトグラフシグナルと全範囲スペクトルを同時に計算し、データシステ ムに送信します。

UV クロマトグラムまたはシグナルは、吸収率データ / 時間軸のプロットで す。波長と帯域幅で定義されます。

- ・ 波長は、検出バンドの中央を示します。
- ・帯域幅は、波長の範囲を定義します。この範囲で吸光度の値の平均が計算され、計測時刻ごとに表示されます。

3 Agilent 1290 Infinity LC システムの最適化 感度を向上させる方法

感度を向上させる方法

例えば、波長が 250 nm で帯域幅が 16 nm のシグナルは、242 nm ~ 258 nm の吸光度データの平均です。また、各シグナルでは、リファレンス 波長とリファレンス帯域幅も定義できます。リファレンス波長の中心にあ るリファレンス帯域幅から計算される平均吸光度が、シグナル波長の相当 する値から減算されて、出力クロマトグラムを生成します。

シグナルの波長と帯域幅は、以下の項目で最適化されるように、選択でき ます。

- ブロードバンドユニバーサル検出
- ナローバンド選択検出
- 特定の検体に対する感度

ブロードバンドまたはユニバーサル検出には、その範囲の吸光度を持つ任 意の試料を検出するための広い帯域幅が必要です。例えば、200 nm ~ 300 nm のすべての吸光分子を検出するには、250 nm のシグナルに帯域幅 100 nm を設定します。欠点は、これらの分子のいずれに対しても感度が最 適化されないということです。ナローバンドまたは選択検出は、最も多く 使われます。特定の分子に対する UV スペクトルが調べられ、適切な極大 波長が選択されます。溶媒が強く吸収する範囲は、可能ならば、回避しま す(メタノールの場合 220 nm 以下、アセトニトリルの場合 210 nm 以下)。例えば、『63 ページ 図 23』では、アニス酸は 252 nm で適切な極大波 長を持ちます。4 nm ~ 12 nm のナロー帯域幅は、一般に、良好な感度を 持ち、ナロー範囲の吸光度に適しています。

ナローバンドは、特定の分子の感度に対して最適化されます。帯域幅が広 がるにつれてシグナルが減衰しますが、同時にノイズも減衰するので、ベ ストな S/N の最適値が存在します。およその目安として、この最適値は、 しばしば、UV スペクトルの半分の高さにある自然な帯域幅に近くなりま す。アニス酸の例では、30 nm です。

分子に対する感度を上げるため、分析波長には、通常、最大波長を設定し ます。検出器は、多くのアプリケーションで、2 AU 以上に線形に上昇しま す。そのため、濃度に対して広い線形範囲が提供されます。高濃度分析の 場合は、濃度の線形範囲は、波長に最小波長のような吸光度の低い波長を 設定したり、通常低い吸光度の値も含んでいる広い帯域幅を採用したりす ることによって拡張できます。定量分析のために最大および最小波長を使 う方法は従来の UV 検出器で使われていましたが、スペクトルの急激なス ロープ部分を回避するために、移動させるグレーティングの機械的な耐性 が必要でした。ダイオードアレイベースの検出器はこのような制限は持っ

感度を向上させる方法

ていませんが、従来の慣習に従って、最大値や最小値が、スペクトル内の 他の値よりも優先して選択されます。

リファレンス帯域幅は、通常、UV スペクトル上で検体の吸光度が 0 の領 域に設定されます。これを、『63 ページ 図 23』のアニス酸のスペクトルに 示します。このスペクトルは、UV 発色団を含む、大部分の小さな分子で一 般的です。ベストの結果を得るには、リファレンスは、シグナル波長にで きるだけ近く、吸光度はゼロの領域上に、ワイドバンドとなるように設定 します。通常、60 nm ~ 100 nm のリファレンス帯域幅が使われます。デ フォルトのリファレンスは、360 nm で、帯域幅 100 nm です。広い帯域幅 が使われるのは、リファレンスシグナル内のノイズが低減できるからです (統計理論によれば、エラー(この場合はノイズ)は、測定値の平方根に 比例して減少します)。リファレンス帯域幅は、スペクトルの吸光度がゼ ロではない部分までは拡張できないことに注意してください。拡張すると、 結果のシグナルは減衰し、感度も不良になるからです。リファレンス波長 を使用すると、温度変化やグラジエント操作が原因の屈折率の変化による、 クロマトグラムのドリフトまたはうねりを減らすことができます。リファ レンスシグナルの効果は、リファレンスシグナルを加えたか加えなかった かだけが異なり、それ以外は同じシグナルを設定することで容易にテスト できます。吸光度が 0 のスペクトル領域がない場合は、リファレンスシグ ナルは無効にしておくことをお勧めします。



Agilent 1290 Infinity LC システムマニュアルおよびクイックリファレンス

3 Agilent 1290 Infinity LC システムの最適化 感度を向上させる方法

スリット幅(G4212A のみ)

分光器への光透過率かつと光学帯域幅は、開口部 可変絞り入射スリット で制御します。スリット幅の初期設定は 4 nm ですが、この設定は、良好 な全方向パフォーマンスを持っているので、大部分のアプリケーションに 適しています。影響を受けるパフォーマンス特性は、感度、スペクトル分 解能、直線性です。分光器へ入る特定の波長を考えると、その光はダイ オードの小さなバンドに対応し、その幅は入射スリットの幅に比例します。 幅が 4 nm のスリットを使って、この状況を説明します。光は 4 nm の帯 域幅を検出した多数のダイオードに対応します。最小光学分解能は 4 nm となり、ダイオードアレイ (またはデジタル)帯域幅には、4 nm 以上を 設定する必要があることになります。最適の感度の場合は、8 nm の設定で ほとんどすべての入力光を受けることができ、ノイズも最小になりますが、 スペクトル分解能は最低になります。これは、UV スペクトルでは通常問題 にはなりません。その自然な帯域幅が、微細構造を持たず、25 nm よりも 大きいからです。8 nm の光学帯域幅は、4 nm のスリットに比べて、線形 範囲が狭くなります。そのため、バリデーション済みのメソッドでは、必 ずバリデーションで使用したスリット幅を使用することが重要になります。 最適なスペクトル分解能のためには、1nm の設定がベストです。この設定 を使えば、ベンゼンのスペクトルに見られるような微細構造が分離できま す(『64ページ図24』を参照)。溶液スペクトル内にこのような微細構造 を表示する化合物は非常にまれです。光度を下げれば、シグナル内にノイ ズが増えます。ノイズレベルは、波長と使用中の移動相溶媒に依存します。



図 24 1 nm および 4 nm のスリット幅の場合のベンゼン(原理)

注入量とサンプル溶解溶媒は、拡散を制御する上で重要です。ピークの高 さが下がる可能性があるので、注入によるピーク分散を避けるために、化

64

感度を向上させる方法

合物はカラムの上部にある部分が分析対象となるよう、注意する必要があ ります。そのためには、サンプルを、移動相よりも低い溶出度の溶媒組成 で溶解する必要があります。カラムで注入量を増やして検体の濃度を高め ることによって、ピーク高さを高くする方法もあります。

『「注入量を増加させる方法」51ページ図』内のコメントを参照してください。

ピーク幅、レスポンスタイム、データ取込レート

検出器内のピーク幅の設定、レスポンスタイム、取込速度は、すべてリン クしています。利用可能な設定を、『66 ページ 表 8』に示します。最適の 感度を得るため、そして分離で得られた分解能を保つためには、これらの 値を正しく設定する必要があります。

検出器は、内部では、クロマトグラムで必要とされる速度よりも速い速度 でデータ数を取込し、それらを処理して、データシステムで表示できるシ グナルを生成しています。処理の一部には、データを適切な取込速度に下 げる処理も含まれ、そうすることによって、クロマトグラフピークの正確 な描画を可能にしています。大部分の分析測定と同様に、読み取られた値 は平均化され、結果内の誤差を低減しています。検出器は、生データ数を グループ化し、電子フィルタリング処理で必要とされるデータ取込レート で、出力シグナルデータを生成します。結果の取込速度が遅すぎる(オー バーフィルタリング)場合は、ピーク高さが低くなり、ピーク間の分解能 が低下します。速すぎる場合は、データにノイズが混じるようになり、狭 いピークを正確にプロファイルできなくなります。

これらのパラメータは、検出器内のピーク幅設定を使えば、クロマトグラム積分結果のピークが広すぎると認識したこと以外の知識なしに、正しく設定できます。ピーク幅設定には、クロマトグラム内の測定値のうち最も狭いピーク幅を設定する必要があります。広すぎる値を設定すると、ピークは高さが低くなり、幅が広がり(そして、おそらく分解能は低下し)ます。狭すぎる値を設定すると、ベースラインノイズを不必要に増やします。 基本的にソフトウェアはこの値を使って、データ取込レートを設定し、最も狭いピークにまたがって十分なデータ数を収集できるようにします。また、ピーク内では15~25ポイントを目標にします。1290 Infinity DAD は、必要に応じて、最大160 Hz で収集できますが、この場合は、わずか0.1 s 幅のピークにまたがって十分なデータ数を収集できます。レスポンスタイム設定は、フィルタリングを設定するためのもう1 つの方法です。レスポンスタイムは秒単位で測定しますが、これは(分単位で測定した)

感度を向上させる方法

ピーク幅の値の約 1/3 になります。この値には、プロットされるシグナル が、入力シグナルのステップの変化に対応する時間を示します。

表 8 ピーク幅 - レスポンスタイム - データ速度

半値幅 [分] ¹	レスポンス [秒]	シグナル取込速度 [Hz]
< 0.0015625	0.015625	160
> 0.0015625	0.03125	160^{2}
> 0.003125	0.0625	80
> 0.00625	0.125	40
> 0.0125	0.25	20
> 0.025	0.5	10
> 0.05	1.0	5
> 0.10	2.0	2.5
> 0.20	4.0	1.25
> 0.40	8.0	0.625
> 0.85	16.0	0.3125

¹ ユーザーインタフェース内の値は、四捨五入されます。

キャリーオーバを最少にする方法

キャリーオーバを最少にする方法

後続するブランク溶媒分析に、直前に分析した注入のピークが観察される 場合に、キャリーオーバが計測されます。キャリーオーバはアクティブな 注入間で持ち越され、誤差のある結果を生成します。キャリーオーバのレ ベルは、直前にアクティブだった注入の領域の割合を使って、ブランク溶 液内のピーク領域としてレポートされます。Agilent 1290 Infinity オー トサンプラは、流路の注意深い設計とサンプル吸収が最小の物質を使用す ることによって、最小のキャリーオーバに対して最適化されています。ト リプル四重極質量分析装置が検出器の場合でも、キャリーオーバ値 0.002 % を達成する必要があります。オートサンプラのオペレーティング 設定を使うと、システム内に粘りつきやすい化合物を含む、任意のアプリ ケーションのキャリーオーバを最小にするための適切なパラメータを設定 できます。

キャリーオーバを最小にするには、オートサンプラの以下の機能を使用します。

- 内部ニードルの洗浄
- 外部ニードルの洗浄
- ニードルシートのバックフラッシュ
- 注入バルブのクリーニング

ニードルの内部を含む流路は、通常のオペレーション中に絶えずフラッシュされているため、大部分の状況ではキャリーオーバを完全に排出します。自動ディレイボリューム削除(ADVR)機能はディレイボリュームを低減しますが、同時にオートサンプラのフラッシングも抑えてしまうので、 キャリーオーバが問題となる対象化合物では使ってはいけません。

ニードルの外部は、特定の場所にある洗浄バイアルを使って洗浄できます。 また、ニードルは、フラッシュポートを使って洗浄できます。トレイ位置 を指定して洗浄バイアルを選択する場合は、セプタムのないバイアルを選 択する必要があります。また、バイアルにはニードルのサンプルを洗浄す るのに適した溶媒が含まれている必要があります。ニードルが下がる際の 汚れが再びニードルが上がる際に汚れが付く可能性があります。ニードル はバイアル中に複数回浸すことができます。これは少量のキャリーオーバ

キャリーオーバを最少にする方法

を取り除くのには有効ですが、ニードルの外部を効率的に洗浄するにはフ ラッシュポートを使う必要があります。

フラッシュポートはニードルシートの上部背面にあり、ペリスタルポンプ によって洗浄溶媒が供給されます。フラッシュポートは、0.68 mlのボ リュームを持ち、ペリスタルポンプは 6 ml/分を供給します。つまり、フ ラッシュポートボリュームには、7 秒で完全に新鮮な溶媒が満たされるこ とになります。フラッシュポートを選択する場合は、新鮮な溶媒でニード ル外部を洗浄する時間を指定できます。キャリーオーバの問題が少ない通 常の状況では、2 秒または 3 秒の短い時間を指定します。完全な洗浄が必 要な場合は、10 ~ 20 秒 を指定します。ニードルシートを汚さないため に、フラッシュポート内のニードルの外部の洗浄は、標準の手順にしてお くことをお勧めします。ニードルシートが汚れた場合は、手操作でフロー 接続を変更してクリーニングすることによる、バックフラッシュが必要で す。これは、フレキシブルキューブモジュールを使って自動化できる操作 の1 つです。

フラッシュポート、溶媒送液ポンプ、接続チューブは、キャリーオーバを 最小にするために定期的にフラッシュする必要があります。たとえば、毎 日、システムを使う前に、フラッシュポンプを3分間、適切な溶媒を使っ てプライミングします。

キャリーオーバを除去するための他の対策が失敗した場合は、インジェク タバルブの内側に対象化合物が残ったままになっている可能性があります。 そのため、キャリーオーバで問題が生ずる場合は、インジェクタバルブに、 バルブ内の流路をクリーニングするための追加の切り替え動作を設定でき ます。問題の化合物が溶出のために高率の有機相を必要とする場合は、最 後のピークを溶出した後に、注入バルブを高率の有機相に切り替えること をお勧めします。また、移動相の初期条件が安定した後は、注入バルブを 再び切り替えておくことをお勧めします。それにより、バルブのロータ シール内のバイパスグルーブにグラジエントの開始条件が設定されること が保証されます。これは、0.5 ml/分以下の流量の場合は、特に重要です。

サンプル用にニードルの外部がフラッシュポンプから供給される水やアル コールでは完全に洗浄できない場合は、適切な溶媒を含む洗浄バイアルを 使用します。インジェクタプログラムを使うと、洗浄のために複数の洗浄 バイアルを使用できます。

カラムの詰まりを回避する方法

カラムの詰まりを回避する方法

HPLC システムを使う場合は、不注意な使用により、カラムまたはシステム チューブ接続が、部分的に、または全体が詰まってしまうことを回避する ために、慎重な配慮が必要です。システムに取り込む物質が原因となる問 題は、通常、以下の作業を行うことで回避できます。

- 溶媒のフィルタリング
- サンプルのフィルタリング
- 移動相の定期的な交換
- システムから緩衝塩をフラッシュする

粒子状物質の不可避的な原因は、システムそれ自体にあります。HPLC シス テムでは、シール材を磨滅させ、その結果生じた粒子状物質がシステム内 のフリットで捕獲されるため、フリットを日常的に交換する必要がありま す。サブ 2 マイクロンの粒子が充填されたカラムでも、充填剤の漏出を避 けるため、細孔径のフリットが必要です。そのため、これらのフリットが、 サンプル、移動相、または機器自体から発生する粒子状物質で詰まってし まうリスクが増大します。

最善の結果を得るには、以下の使用ガイドラインに従ってください。

- 1 カラムの取り付けと分析は、必ず、カラム上にマークされているフロー 方向を守って行ってください。
- 2 溶媒は、高品質で、クロマトグラフィグレードのものだけを使ってください。
- 3 すべての水性緩衝液とすべてのサンプルは、使用する前に、0.2 µm フィ ルタを使って、フィルタリングを行ってください。
- 4 移動相緩衝液のボトルは、毎日または2日ごとに交換してください。移動相をボトルに継ぎ足さないでください。必ず、新しいボトルを使ってください。
- 5 高緩衝塩移動相(> 50 mM)は、アセトニトリルと組み合わせて使わない でください。沈殿物が生ずる可能性があるからです。
- 6 高濃度緩衝移動相では、シール洗浄オプションを使ってください。

カラムの詰まりを回避する方法

- 7 粒子状物質を捕獲し、カラムの寿命を伸ばすために、インラインフィル タ(1290 Infinity インラインフィルタ、内径 2 mm、部品番号: 5067-4638)を使うことをお勧めします。
- 8 圧力が 10 % 増加したら、フィルタを交換してください。
- 9 カラムを機器に取り付ける前に、移動相を含む緩衝液のポンプ(カラム までの接続部分)をパージし、5 mlの溶媒を使ってフラッシュしてく ださい。
- 10 カラムを互換性のある移動相でフラッシュします。0.1 ml/分(内径 2.1 mm のカラムの場合)、0.2 ml/分(内径 3.0 mm のカラムの場合)、0.4 ml/分(内径が 4.6 mm のカラムの場合)の流量でゆっくりと開始してください。流量を 5 分以上、必要な流量に増やしてください。
- 11 圧力が安定したら、カラムを検出器に接続してください。
- 12 カラムと検出器は、使用する前に、移動相をカラムの 10 倍以上のボ リュームを使って平衡を保ってください(カラムボリュームについて は、『44 ページ 表 7』を参照)。
- 13 過圧力は避けてください。シーケンスを開始する前に、グラジエントの 圧力範囲をチェックしてください。
- 14 溶解したサンプルの有機含有量が多い場合は、注入プログラムを使っ て、サンプルの前後に弱い溶媒のプラグを追加し、高濃度緩衝液移動相 内の沈殿のリスクを低減することを考えてください。



4

Agilent 1290 Infinity LC システムマニュアルおよびクイック リファレンス

. システムの設定とインストール

ソフトウェアのインストール 72
モジュールの設置 74
スタック コンフィグレーションの最適化(バイナリ LC シ ステム) 74
スタックコンフィグレーションの最適化(クォータナリ LC システム) 79
ポンプのプライミング 84
ポンプのパージ 87
モジュール間のフロー接続 90
ネットワークへの統合 90

この章では、ソフトウェアのインストール、スタック構成、システ ム運用の準備について説明します。



ソフトウェアのインストール

ソフトウェアコントローラとデータシステムのインストー ル

ソフトウェアのインストール手順については、1290 Infinity Diode Array Detector Manual とソフトウェアマニュアルを参照してください。

Agilent Lab Advisor ソフトウェアのインストール

Agilent Lab Advisor ソフトウェアのインストールについては、Lab Advisor DVD のソフトウェアマニュアルを参照してください。

Agilent Lab Advisor は、従来は ChemStation ソフトウェアに含まれてい た診断機能に代わるものです。

Agilent Lab Advisor は、ラボ内の機器をリアルタイムで常時監視するための Windows® ベースのアプリケーションであり、高度なカウンタを使って保守とサービスの必要性を自動的に通知することで生産性を向上させます。この機能を使うと、問題点は結果に影響を与える前に修復できます。このソフトウェアには、多くのユーザー情報とマニュアル、一連のカルキュレータ、機器の設定、キャリブレーション、保守を支援するツール、正常なパフォーマンスを検証するためのテストおよび診断ルーチンが含まれています。また、Agilent Lab Advisor は、発生する可能性のある機器障害について、フィードバックと解決策を備えています。このソフトウェアは、Agilent データシステムの有無に関わらず使用できます。

このソフトウェアの監視対象は、以下のとおりです。

- LC モジュールのステータス
- 早期保守フィードバック(アップグレードまたは交換の必要性を判定するため)

また、以下の機能も持っています。

• 診断テストの自動化
ソフトウェアのインストール

- 電源が投入され、PC またはラボのネットワークに接続されている、サポート対象の LAN ベース機器の識別
- 一般的な機器不具合について、部品交換やトラブルシューティング作業 を自動的に診断

4 システムの設定とインストール モジュールの設置

モジュールの設置

モジュールの設置方法については、個々のモジュールのマニュアルを参照 してください。これらのマニュアルには、仕様、保守、部品についての説 明も記載されています。

スタック コンフィグレーションの最適化 (バイナリ LC シ ステム)

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(『75ページ 図 25』と『76ページ 図 26』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。

システムの設定とインストール 4 モジュールの設置



図 25 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)

4 システムの設定とインストール モジュールの設置



Agilent 1290 Infinity LC システムマニュアルおよびクイックリファレンス

システムの設定とインストール 4 モジュールの設置

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(『77 ページ 図 27』および 『78 ページ 図 28』を参照)。



図 27 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)

Agilent 1290 Infinity LC システムマニュアルおよびクイックリファレンス

4 システムの設定とインストール モジュールの設置



AC 電源

図 28 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)

システムの設定とインストール 4 モジュールの設置

スタックコンフィグレーションの最適化(クォータナリ LC システム)

1スタック構成

以下のコンフィグレーションで Agilent 1290 Infinity クォータナリ LC システムのモジュールをインストールし、確実に最適な性能が得られるよ うにしてください(『80ページ 図 29』と『81ページ 図 30』を参照)。こ のコンフィグレーションでは、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を最小限に抑えます。

Agilent 1290 Infinity クォータナリポンプは、常にスタックの最下部に 設置する必要があります。 4 システムの設定とインストール モジュールの設置



クォータナリ ポンプを装備した 1290 Infinity の推奨スタック構 図 29 成(前面図)

システムの設定とインストール 4 モジュールの設置



図 30 クォータナリ ポンプを装備した 1290 Infinity の推奨スタック構 成(背面図)

Agilent 1290 Infinity LC システムマニュアルおよびクイックリファレンス

4 システムの設定とインストール モジュールの設置

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、2 スタッ ク コンフィグレーションを推奨します。重いモジュール 両方 (1290 Infinity ポンプとサーモスタット)を各スタックの最下部に設置し、ス タックが高くなるのを避けられます。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。(『82 ページ 図 31』と『83 ページ 図 32』を参照)。



図 31 クォータナリ ポンプを装備した 1290 Infinity の推奨 2 スタック 構成(前面図)

Agilent 1290 Infinity LC システムマニュアルおよびクイックリファレンス

システムの設定とインストール 4 モジュールの設置



AC 電源

図 32 クォータナリ ポンプを装備した 1290 Infinity の推奨 2 スタック 構成(背面図)

4 システムの設定とインストール モジュールの設置

ポンプのプライミング

この手順が必要になるのは、以下の場合です。

- ポンプを初めて使う
- 1 つ以上のインレットチューブが気泡を含んでいる、または何らかの理 由で乾いている

プライミングの目的は、ポンプとインレットチューブからすべての気泡を 取り除くことです。

2-propanol (別名: イソプロパノール、イソプロピルアルコール、IPA) は、乾いているポンプのプライミングに最適な溶媒です。大部分の順相ま たは逆相の溶媒と親和性があるという利点を持ちます(不明な場合は、溶 媒混和性の表を参照)。

有毒、可燃性および有害な溶媒、サンプル、試薬 警告

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。
- 溶媒の容器を空にして運用することを避けるため、Bottle Filling モニタを 注記 使います。シリンジを使って、溶媒チューブを満たします。ポンプは空気は排 出しません。
- 手順を開始する前に、すべての溶媒ラインが、ポンプ設置マニュアルの記載に 注記 従って、ポンプに接続されていること、および廃液チューブが溶媒廃液コンテ ナに取り付けられていることを確認します。ポンプがコンピュータソフトウェ アまたは Instant Pilot コントローラで制御されていること、および流量に ゼロが設定されていることを確認します。

- モジュールの設置
- 1 チャンネル A とポンプヘッド A をプライミングするために準備するに は、以下の手順を実行します。
 - a ポンプをプライミングするため、溶媒の各容器を約 150 ml の HPLC グレードの 2-propanol で部分的に満たし、溶媒チューブのガラス フィルタの端を容器に取り付けます。
 - b ポンプヘッド A のインレットチェックバルブに接続しているチューブを取り外します。これは真空デガッサチャンネル A のアウトレットチューブです。
 - **c** ねじ込み式アダプタが付いたプライミングシリンジを、チューブに接 続します。
 - d すべての気泡が取り除かれるまで、溶媒をゆっくりとチューブから吸 引します。
 - e 直後に、チューブからシリンジとアダプタを取り外し、ポンプヘッド A のインレットチェックバルブにそのチューブを接続します。
- 2 溶媒チャンネル B とポンプヘッド B に対して、上記の手順を繰り返し ます。
- 3 ポンプに溶媒切り替えバルブ(SSV)が取り付けられている場合は、残り の空の溶媒チューブの各々に対して、以下の手順を実行します。
 - a 溶媒切り替えバルブから空のチューブを取り外します。ねじ込み式ア ダプタが付いたプライミングシリンジを、チューブに接続します。
 - b すべての気泡が取り除かれるまで、溶媒をゆっくりとチューブから吸 引します。
 - c 直後に、チューブからシリンジとアダプタを取り外し、溶媒切り替え バルブにそのチューブを再接続します。
- 4 ChemStation ソフトウェアで、システムダイアグラムのポンプセクションを右クリックして、コンテキストメニューから プライミング オン を 選択します (パージオン コマンドと間違えないでください)。他のコントローラにも同様の機能があります。

この段階で、パージバルブは流路 A と B を廃液ポジションに切り換え、 それと同時に、両方のチャンネルをプライミングします。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。これが 20 回繰り返された後、 プライミングプロセスは完了します。パージバルブは、流路を切り換えて システムに戻します。

モジュールの設置

この段階で各チャンネルを 30 ml の 2-propanol を使ってパージすること をお勧めします。『「ポンプのパージ」87ページ 図』の手順に従います。

ポンプのパージ

警告

1290 Infinity バイナリポンプのパージ手順について説明します。1290 Infininty クォータナリポンプの場合も、同じ方法で実施できます。

- ポンプが初めてプライミングされた後。
- システムを使用する前にポンプを新しい溶媒でパージするとき、または 溶媒を交換するとき。
- ポンプが数時間以上アイドルになっていたとき(溶媒ライン内に空気が 拡散している可能性があるので、パージすることをお勧めします)。
- 溶媒ボトルが満たされ、システムに新しい溶媒を満たすためにポンプで パージが必要になったとき。異なる溶媒を使う場合は、新しい溶媒が前 の溶媒と混和性があることを確認してください。必要に応じて、coと 混和性がある溶媒(多くの場合に、イソプロパノールが適しています。 溶媒混和性の表をチェックしてください)が使われる中間ステップを 使ってください。

ポンプへ至るインレットチューブは、すでに溶媒で満たされている必要が あります。インレットチューブが部分的または全体的に乾いている場合は、 プライミング手順の全体を実行します(『「ポンプのプライミング」84 ページ 図』を参照)。

パージバルブを使うと、両方のポンプヘッドを同時に廃液系に接続し、そ れぞれの最大流量 5 ml/min でパージできます。50/50 混合に対する合計 パージ流量は 10 ml/min となります。

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。

モジュールの設置

 パージバルブを制御するための設定ページにアクセスするには、ポンプ セクションを右クリックして、コンテキストメニューから コントロー ル を選択します。

あるいは、Instrument > More 1290 Infinity BinPump > Control を選 択します。

Control	×
Pump	
	 Off Standby
Seal Wash	
 Off Single Wash Periodic 	Duration 1.0 0 min Period 2.0 0 min
	on for 0.2 timin
Automatic Turn On	
🔲 Tum on at	Monday, July 06, 2009 11:00:00 🗘 👻
Purge	
© On ◉ Off	Duration 1.00 + min. Flow 0.000 + mL/min.
Prime © On Off	
	<u>O</u> k <u>C</u> ancel Help

- 2 パージ セクションで、以下のパラメータを設定します。
 - 時間:6 分
 - 流量: 10 m1/分
 - 混合 B: 50 %

モジュールの設置

混合 A は、自動的に 50 % が仮定されます。On/Off ボタンは、オフ の 設定のままにしておきます。OK をクリックして終了します。

ポンプセクションを右クリックして、コンテキストメニューからパー
 ジオンを選択します。

| **パージ オン** を次の項目 **プライミング オン** と間違えないでください。

注記

この段階で、パージバルブは流路をパージポジションに切り換え、両方の チャンネルを 5 ml/min の流量で廃液系に同時にパージします。それぞれ の所要時間は 6 分です。設定した所要時間が経過すれば、パージ流量はオ フとなり、バルブは流路をシステムに戻します。メソッドで設定されてい る流量と混合が再開します。この例題では、メソッドのフローはゼロに設 定されたままです。手順 1~2 の **コントロール** ページのパージ設定が維 持されるため、再びパージプロセスが必要になった場合は手順 3 から開始 できます。

ポンプが最初からプライミング済みで、propan-2-o1 を使ってパージされ ていた場合は、溶媒を水やメタノールのような移動相溶媒と交換できます。 溶媒を交換した場合は、パージ手順を繰り返します。パージが終了した直 後は、ポンプ内の溶媒は脱気されていません。したがって、少なくとも 10 min間はシステムの運転を続けて、溶媒を脱気する必要があります。

モジュールの設置

モジュール間のフロー接続

モジュールを接続する場合は、各キャピラリと溶媒を含むカラムを、流路 内の次のコンポーネントに接続する前に、フラッシュする必要があります。

- Jet Weaver ミキサーのアウトレットを、内径 0.12 mm のステンレス製 フレキシブルキャピラリ(カラーコードラベルは赤)を使って、オート サンプラに接続します。これは、オートサンプラのインジェクションバ ルブのポート 1 に接続する必要があります。
- 2 内径 0.12 mm のステンレス製のフレキシブルキャピラリを、オートサン プラのインジェクションバルブのポート 6 から、温度調節機能付きカラ ムコンパートメントに接続します。キャピラリは、ディレイボリューム が最小になるように、低拡散ボリューム熱交換器に直接接続します。切 り替えバルブが取り付けられている場合は、それに接続します。
- 3 カラムのアウトレットを、1290 Infinity 検出器内の Max-Light カート リッジセルのインレット(CELL-IN とマークされている、左側の接続) に接続します。
- 4 廃液チューブを、1290 Infinity 検出器内の Max-Light カートリッジセルのアウトレット(CELL-OUT とマークされている、右側の接続)に接続し、アウトレットを廃液収集コンテナ内に配置します。

ネットワークへの統合

システムをネットワークに接続する方法については、モジュールのユー ザーマニュアルを参照してください (LAN Configuration の章)。



5

Agilent 1290 Infinity LC システムマニュアルおよびクイック リファレンス

クイックスタートガイド

クイックスタートガイドについて 92 システムの準備 93 システムをオンにする 93 デフォルトメソッドの読み込み 94 オンラインプロットの構成 95 ポンプのパージ 97

Method and Run Control (メソッド & ランコントロール) ビュー
内のデータ取得 98
テスト用の混合物および ZORBAX RRHD カラム用のメソッドパ
ラメータ 98
メソッドの設定 100
シングルラン用のメソッドの実行 103
メソッドの高速実行 105

データ分析 107
 データ分析ビュー 108
 シグナルの積分 109
 レポートの指定 111

この章では、1290 Infinity LC システムを使ったデータ取込とデー タ分析について説明します。



クイックスタートガイドについて

クイックスタートガイドについて

この章では、Agilent 1290 Infinity LC システムを運用する方法について 説明します。この章は、インストール後に初めて分析を行うためのクイッ クスタートガイドです。チュートリアルの例題として使ったり、システム 全体の機能を確認するために使います。また、メソッド パラメータの詳し い説明も含んでいます。

この例題では、Agilent 1290 Infinity LC システムに用意されているカラ ムと点検サンプルを使って、設定と分析を行う方法について説明します。 この例題では OpenLAB CDS ChemStation エディションのメニューとコマン ドを使いますが、同等の機能は、OpenLAB CDS EZChrom エディション、イ ンスタントパイロットコントローラ、MassHunter ソフトウェアなどの代替 制御オプションでも用意されています。

注記 ここでは、システムはインストール済みで動作中であり、初期プライミングも 終わっていると仮定します(『「ポンプのプライミング」84ページ 図』を参照)。UV ランプは、少なくとも、定量解析を行う 30 分前にオンにしておく必要 があります。

システムの準備

システムをオンにする

システムの準備が完了していない場合、つまりレディ状態ではないソフト ウェアがある場合は、以下の手順を実行します。

- コンピュータシステムをオンにして、Windows デスクトップが表示されるのを待ちます。
- 各 LC モジュールの左下にあるボタンを使って、電源をオンにします。
 ボタンの中央には、緑のランプが見えます。
- 3 アイコンをクリックして、コントロールソフトウェアを起動します(構成済みの場合)。あるいは、スタート > すべてのプログラム > Agilent Technologies > OpenLab > OpenLab コントロールパネル を選択します。 機器の下のナビゲーションウィンドウで該当する機器を選択し、オン ライン起動をクリックします。

メソッド & ランコントロール ビューに、ChemStation ソフトウェアが 表示されます。モジュールは、最初はスタンバイモードでノットレディ 状態です。ただし、直ちに初期化が行われ、レディ状態となるオートサ ンプラだけは例外です。

4 各モジュールを個別にオンにするには、対応するアイコンを右クリック して、コンテキストメニューから[モジュール名]に切り替えオン を 選択します。

あるいは、システムダイアグラムの右下にある システム オン/オフ ボ タンをクリックして、システム内のすべてのモジュールを一斉にオンに します。時間が経過して設定ポイントに到達すると、システムステータ スは、Not Ready(黄色)から Ready(緑色)に変化します。

デフォルトメソッドの読み込み

ChemStation には、デフォルトのメソッド **DEF_LC.M** があります。このメ ソッドは、最初の実行時、または新しいブランクテンプレートが必要に なった時点で、読み込まれます。このメソッドには、すべてのモジュール のデフォルト設定が含まれています。

この手順では、メソッド DEF_LC.M を読み込みます。このメソッドを使って、すべてのパラメータにデフォルト設定を設定します。またはブランクのメソッドテンプレートを取得して、新しいメソッドを設定します。

- **1** ChemStation の メソッド & ランコントロール ビューに移動します。
- 2 メニューバーで メソッド > 新しいメソッド... を選択し、コンテキス トメニューから DEF_LC.M を選択します。

あるいは、メニューバーの下の メソッド読み込み アイコン 崎 を使ったり、ナビゲーションペイン内の メソッド タブのメソッド名 DEF_LC. M をダブルクリックします。



デフォルトのメソッド (DEF_LC. M) は、一連のデフォルトパラメータを 持っており、これらを変更することによって新しいメソッドを作成します。 たとえば、流量にゼロを設定し、メソッド情報 と メソッド履歴 をブラン クにします。

注記

このメソッドには、新しいパラメータは上書きできないことに注意してください。したがって、上書き保存 をクリックすれば、名前を付けて保存 機能にリダイレクトされるので、異なるメソッド名を入力する必要があります。

オンラインプロットの構成

 オンライン プロット ウィンドウが表示されない場合は、以下の手順を 実行します。View > Online Signals > Signal Window 1 をクリックす ると、ウィンドウが表示されます。

	Online Plot						
	DAD1A, DA	D: Signal A					
	mAU - 60 - 40 - 20 -						
	0	0.5	1	1.5	2	2.5	min
÷		Change			Adjust		4

2 オンライン プロット ウィンドウ内に必要なシグナルを構成するには、 変更… をクリックします。

Edit Signal Plot	
<u>A</u> vailable Signals	Selected Signals
PMP1B, BinPump: Pressure PMP1C, BinPump: Flow PMP1D, BinPump: Solvent Ratio A PMP1E, BinPump: Solvent Ratio B	DAD1A, DAD: Signal A
THMIA, TCC: Left Temperature THM1B, TCC: Right Temperature DAD1B, DAD: Signal B DAD1C, DAD: Signal C	<- <u>R</u> emove
Window <u>×</u> -axis 3 min	DAD1A, DAD: Signal A Type: acquired ⊻-axis range: 100 ★ mAU
🗖 draw zero line	☐ auto y-adjust Offset: 10 📕 %
Fraction Collector	Method Settings Apply to Method
ок	Cancel Help

シグナルプロットの編集 設定ページが表示されます。



- 3 使用可能シグナル ボックスで必要なシグナルを強調表示し、追加 をク リックして、それらを **シグナル選択** ボックスに移します。
- 4 各シグナルに対して個々の設定を行うには、シグナル選択 ボックス内 のシグナルを強調表示し、ページの下半分で必要な値を設定します。

検出器シグナルの他に、温度や圧力のようなパラメータのトレース結果もプ 注記 ロットできます。Apply to Method を使って、このページ内の設定をメソッド に保存します。

> オンラインプロット ウィンドウは、電子グラフ用紙のような働きをしま す。つまり、検出器の出力その他の出力パラメータを継続して記録します。 シグナルはウィンドウの右から左に描画されます。最高で 90 分 の過去の データにアクセスできます。これはベースラインをチェックし、前の注入 を観察するのに役立ちます。X 軸と Y 軸のスケールは、各軸上の上 / 下ボ タンを使って直接調整できます。

> オンライン プロット ウィンドウの 調整 ボタンを使えば、 選択したシグ ナルの現在のポイントをゼロラインに移動できます。選択したシグナルは、 Y 軸のラベルの色で示されます。シグナルの上、またはプロット上部の関 i 車シグナルの説明をクリックすることによって、特定のシグナルを選択で きます。

Balance ボタンをクリックすると、検出器の出力信号がゼロになります。

オンライン プロット ページを変更しても、個々のデータファイルに保存され 注記 ているデータには一切影響はありません。

ポンプのパージ

以下の状況の場合は、ポンプをパージします。

- ポンプが初めてプライミングされた。
- ポンプを、システムを使う前に、新しい溶媒でパージするとき、または 溶媒を交換するとき。
- ポンプが数時間以上アイドルになっていたとき(溶媒ライン内に空気が 拡散している可能性があるので、パージすることをお勧めします)。
- 溶媒の容器が満たされ、システムに新しい溶媒を満たすためにポンプで パージが必要になったとき。異なる溶媒を使う場合は、新しい溶媒が前 の溶媒と混和性があることを確認してください。必要に応じて、2 つの 溶媒と親和性がある溶媒(多くの場合に、イソプロパノールが適してい ます。溶媒混和性の表をチェックしてください)が使われる中間ステッ プを使ってください。

パージ手順についての詳細は、『「ポンプのパージ」87ページ 図』を参照 してください。

Method and Run Control (メソッド & ランコントロール) ビュー内のデータ取得

Method and Run Control (メソッド & ランコントロール) ビュー内のデータ取得

1290 Infinity バイナリポンプのすべてのメソッドの手順と設定について 説明します。1290 Infininty クォータナリポンプの場合も、同じ方法で実 施できます。

テスト用の混合物および ZORBAX RRHD カラム用のメソッド パラメータ

1290 Infinity LC システムには、ZORBAX RRHD Eclipse Plus C18 (1.8 µm、2.1 mm x 50 mm) カラムと、ここでの例の手順で使用するための テスト用の、phenone を含む、混合物が付属しています。

Phenones Test Mix (部品番号 5188-6529) は、各々が水 / アセトニトリ ル (65/35) に溶解する 100 ng/µl のコンポーネントを 9 つ持っていま す。9 つのコンポーネントは以下のとおりです。

- アセトアニリド
- アセトフェノン
- プロピオフェノン
- ・ ブチロフェノン
- ・ ベンゾフェノン
- ・ バレロフェノン
- ヘキサノフェノン
- ヘプタノフェノン
- オクタノフェノン

このテスト用混合物の分離用のメソッドパラメータを、『99 ページ 表 9』 に示します。

Method and Run Control (メソッド & ランコントロール)ビュー内のデータ取得

モジュール	パラメータ	設定
ポンプ	溶媒 A	水
	溶媒 B	アセトニトリル
	流量	0.4 m1/分
	最初の混合	60 % A, 40 % B
	グラジエントタイム テーブル	4 分後に 20 % A、80 % B
	ストップタイム	5 分
オートサンプラ	注入量	1 µl
	ニードル洗浄	フラッシュポート、 6 秒間
TCC	カラム	ZORBAX Eclipse Plus C18 1.8 µm, 2.1 mm x 50 mm 内径
	温度	40 °C
検出器	シグナル A	250 nm、バンド幅 100 nm、リファレンス 360 nm、バンド幅 100 nm
	ピーク幅	0.025 分(10 Hz)
	スペクトル保存	すべて

表 9 最初の分離分析用のメソッドパラメータ

Method and Run Control (メソッド & ランコントロール) ビュー内のデータ取得

メソッドの設定

この節では、テスト用混合物の条件を用いて分析用のメソッドの条件を短時間で設定する方法を示します。利用可能なすべてのパラメータについての詳細は、『「メソッド全体の編集を使ったメソッドの設定」119ページ図』の付録を参照してください。

デフォルトのメソッド DEF_LC.M は、新しいメソッドを準備するためにすで に読み込み済みです。この段階で主要なパラメータを編集して、新しいメ ソッドを作成できます。ここで説明する例題の分離に対して、『99ページ 表 9』にリストした条件を設定します。

各モジュールの メソッド ページにクイックアクセスするには、モジュールのシステムダイアグラム内を右クリックして、コンテキストメニューから メソッド...を選択します。

各モジュールをこの方法で設定します。

- 2 ポンプ領域を右クリックして、コンテキストメニューから メソッド... を選択します。
 - a 1290 Infinity Binary Pump の メソッド ページに、以下のパラメー タを入力します。
 - 流量: 0.4 m1/分
 - 溶媒 A: 圧縮率のドロップダウンリストから、水 を選択します。
 - 溶媒 B: チェックボックスを選択し、溶媒 B をアクティブにします。
 - %B: 初期值 40 %
 - ストップタイム:5分
 - 最大圧力リミット値: 1200 bar
 - **b** + 符号をクリックして、**タイムテーブル** を表示します。
 - c ラインを追加し、溶媒混合の変更 を選択して、%B に 80 % を設定し ます。
 - d 他のパラメータは、デフォルトの設定のままにします。OK をクリックしてウィンドウを終了します。
 変更がポンプモジュールに送信されます。

Method and Run Control (メソッド & ランコントロール) ビュー内のデータ取得

- 3 オートサンプラ領域を右クリックして、コンテキストメニューから メ ソッド・・・ を選択します。
 - a 1290 Infinity オートサンプラ の メソッド ページに、以下のパラ メータを入力します。
 - 注入量: 1.0 µl
 - ニードル洗浄を用いた注入
 - モードフラッシュポート、時間:6 秒間
 - b 他のパラメータは、デフォルトの設定のままにします。OK をクリックしてウィンドウを終了します。
 変更がオートサンプラモジュールに送信されます。
- 4 温度調節機能付きカラムコンパートメント(TCC)領域を右クリックして、コンテキストメニューからメソッド...を選択します。
 - a 1290 Infinity TCC の メソッド ページに、以下のパラメータを入力 します。
 - 左側温度 40 ℃
 - 右側温度は左側と同じに
 - b 他のパラメータは、デフォルトの設定のままにします。OK をクリッ クしてウィンドウを終了します。

変更が TCC モジュールに送信されます。

- 5 ダイオードアレイ検出器領域を右クリックして、コンテキストメニュー から メソッド... を選択します。
 - a 1290 Infinity DAD の メソッド ページに、以下のパラメータを入力 します。
 - 使用シグナル:チェックボックスのチェックを外して、シグナルA
 以外のすべてのシグナルをオフにします。
 - シグナル A: 250 nm、バンド幅 100 nm、リファレンス 360 nm、バ ンド幅 100 nm
 - ピーク幅: 0.012 分 (0.25 秒間 レスポンス、20 Hz)
 - b 詳細設定 セクションで、スペクトル保存 に 全て を設定します。
 - **c** 他のパラメータは、デフォルトの設定のままにします。**OK** をクリッ クしてウィンドウを終了します。

Method and Run Control (メソッド & ランコントロール) ビュー内のデータ取得

変更が DAD モジュールに送信されます。

6 この段階で、必要なモジュールパラメータのすべてに値が設定されました。Method > Save Method As... を選択して、メソッドを新しい名前で保存します。

ChemStation ではメソッドを $DEF_LC.M$ として保存することは許されて いません。したがって、デフォルトのメソッドテンプレートが変更され ることはありません。

7 システムが平衡化するまで、少なくとも、10分待ちます。そして、分析を開始する前に、オンラインプロット内のベースラインが安定化していることを確認します。

Method and Run Control (メソッド & ランコントロール) ビュー内のデータ取得

シングルラン用のメソッドの実行

この節では、前の項で入力した条件を使って、テスト用混合物のシングル ランを実行する方法を示します。

ChemStation での分析は、次の 2 つのモードで実行できます。

- ランメソッド たとえば、メソッド開発中のシングルラン。現在設定 されているパラメータ設定を使用する。
- シーケンス開始 複数のバイアルからの一連の注入を自動化。複数のメ ソッドを使う場合もある。詳細は、ChemStation のマニュアルを参照し てください。
- 1 ラン メソッド タスク選択 アイコン 🧧 をクリックします。
- 2 必要なメソッド条件が現在読み込まれていない場合は、メニューバーの 下にある メソッド > メソッド読み込み またはアイコン [№] を選択し、 それらを読み込みます。
- 注記 読み込んだメソッドを変更してまだ保存していない場合は、メソッドステータ スアイコンに黄色のアスタリスクが付けられて、その状況が通知されます。変 更されたパラメータを保存しなくても、注入は行えます。ChemStation は、取 込パラメータのコピーをデータファイル ACQ.TXT に必ず保存します。そのた め、オリジナルの取込パラメータが確実に維持されます。
 - **3** サンプルバイアルをポジション 1 に置きます。ここが、サンプルトレイ の右側の 10 x 2 ml のバイアルポジションの前面ポジションです。
 - 4 ランコントロール > サンプル情報 を選択し、サブディレクトリ 名 (オ プション)、ファイル名、サンプルの ロケーション (バイアル 1)、サンプル名、そして任意の コメント を入力します。
 - 5 システムの平衡化が完了していてベースラインが安定化していれば、サンプル情報ページ内のランメソッドをクリックして、注入を開始します。または、OKをクリックし、準備が完了すれば、システムダイアグラムの上にあるシングルラン開始ボタンをクリックします。
 - 6 注入が行われて、オンライン プロット にクロマトグラムが表示されま す。ストップタイム に到達すると、データ取得が停止します。

Method and Run Control (メソッド & ランコントロール) ビュー内のデータ取得

クロマトグラムは『105ページ 図 33』のように表示されます。ここで は時間軸が短くなっていますが、このクロマトグラムが次項で説明する 4 倍速い条件の下での分析だからです。

Method and Run Control (メソッド & ランコントロール) ビュー内のデータ取得

メソッドの高速実行

最初の練習は、標準システムで達成できる圧力下で行いました。次に、流 量を増加させ、グラジエントを高速分離用に調整します。

- 1 前の項と同様な方法でメソッド条件を編集し、次のように変更します。
 - 流量: 1.6 ml/分
 - グラジエント:グラジエントは、最初の分析と比較してグラジエント スロープのボリュームが変わらないように変更します。流量を4倍に 増やしたので、グラジエント時間も4分の1にします。すなわち、グ ラジエント時間に、1分を設定します。
 - ストップタイム: 1.25 分
- 2 メソッドを新しい名前で保存します。
- 3 サンプル情報 で新しいファイル名が使われていることを確認します。
- 4 ベースラインが安定化し平衡化すれば、シングル サンプル開始 ボタン をクリックして分析を開始します。

クロマトグラムは下図のようになります。実行時間は、約1分です。



題のクロマトグラム

Method and Run Control (メソッド & ランコントロール) ビュー内のデータ取得

この分離は実際にはこれらの条件用に最適化されていません。メソッドを 最適化すれば、システムの実行をさらに高速にすることができます。高速 化に貢献している項目は、以下のとおりです。

- サンプルの濃度を希釈率 1:10 で薄めています。
- グラジエントの範囲を増やしています。
- 温度を上げています。
- ピークのスペクトルを調べて、適切なナローバンド検出機能を選択しています。

最適化についての詳細は、『「Agilent 1290 Infinity LC システムの最適 化」39ページ 図』を参照してください。

データ分析

ChemStation 内のメソッドには、データ取得(システムの制御)用とデー タ分析(定性的定量的結果を取得するためのデータ処理)用のすべてのパ ラメータが含まれています。この項では、データ分析の積分とレポートを 概観し、この章の前半で生成された分離が積分されて出力される状況につ いて説明します。定量分析のためのキャリブレーションの使用を含む、 データ分析についての詳細は、ChemStationのマニュアルを参照してくだ さい。



図 34 データ分析ビュー

データ分析ビュー

データ分析 ビューにクロマトグラムを表示するには、以下の手順を実行します。

- 1 ChemStation をオフライン起動します。
- 画面の左下にある[データ分析]をクリックします(『107ページ図34』 を参照)。
- 3 ナビゲーションパネルで、データファイルを含むデータディレクトリを 探します。すべての単一注入が、シングル ラン のサブセットとして表 示されます。シングル ラン をダブルクリックし、これらのデータファ イルをナビゲーションテーブルに読み込みます。
- 4 ナビゲーションテーブル内のファイルを選択し、それをダブルクリックして、クロマトグラムをビューアに読み込みます。
クイックスタートガイド 5 データ分析

シグナルの積分

1 [積分タスクツール]を選択します(下図参照)。下図では、積分 アイコンと 積分イベント テーブル設定 アイコンが強調表示されています。

ngle Runs: PHENONES								
📙 Integration 🎂 Calibration 🗔 Signal 🔣 Purify	/ 🍐 Spectrum							
📠 🕅 Report: Short 📄 🗋 🖉	i 🗇 🚯							
DAD1 A, Sig=242ONE-MIX-010.D) 💌 🛃 🕀 😪 📘) 🕓 🚑 🖄 ∫	🔊 🔺 📉	<u>ю</u> ,	-				
◈觱із=⋺•Щ┇⊘	DAD1 A, Si	g=242,4 Ref=360,10	00 (PHENONES\P	HENONE-MIX-0	010.D)			
Method Manual Events Final Events For All Signals:	mAU 300 260	2 <u>1</u>	0.192	0.377	0.422	0.487	10918	0.617
Integration Events Value Targent Skim Mode Standard Tail Peak Skim Height Ratio 0.00 Front Peak Skim Height Ratio 0.00 Skim Valley Ratio 20.00 Baseline Concetion Classical Peak to Valley Ratio 500.00	200 - 160 - 100 -			0348				
Specific Events For Signal:			0.2		0.4		0.6	
DAD Default								
-	[•]							
	# Time	Area	Height	Width	Area%	Symmetry		
Time Integration Events Value	1 0.122	281.7	208.1	0.0208	10.633	0.518		
Initial Slope Sensitivity 5	2 0.192	292.2	229.5	0.0189	11.030	0.509		
Initial Peak Width 0.05	3 0.279	257.5	239.1	0.0166	9.720	0.545		
Initial Area Reject 5	4 0.349	24.6	16.8	0.0201	0.929	2.05		
Initial Height Reject 1	5 0.377	232.4	261	0.0144	8.771	0.702		
Initial Shoulders OFF	6 0.422	239	296	0.0124	9.023	0.661		
	7 0.487	257.8	295.9	0.0132	9.730	0.749		
	8 0.601	337.3	337.4	0.0147	12.731	1.421		

- 2 積分イベント テーブル設定 アイコンをクリックして、テーブルを下図 のように開きます。
- 3 グラジエント分析用に、ベースライン補正 に 詳細設定 を設定します。
- 4 スロープ感度 に 50 を設定します。数字を大きくすれば、急なピークを 積分し、緩やかなピークを無視します。
- 5 ピーク幅の値に、対象のピークの中で最も狭いピークを設定します。この例では、約 0.01 です。
- 6 最小のピークをリジェクトするために、面積リジェクトと高さリジェク トを設定できます。
- **7 積分** アイコンをクリックし、これらの新しい設定を用いて、結果を アップデートします。



8 緑のチェックマークアイコンを使って、イベント テーブルを終了しま す(下図参照)。



クイックスタートガイド 5 データ分析

レポートの指定

1 メニューバーで レポート > レポート条件 をクリックすると、下図に示 すウィンドウが表示されます。

Specify Report: Instrument 1	
Reporting settings Quantitation settings	
Report mode Use Classic Reporting	
Style Report style: Short Quantitative result sorted by: Signal Sample info on each page Add fraction to Add chromatogram output Add summed Add sample custom fields to sample info Add compount Report layout for uncalibrated peaks O to not	Chromatogram output Pottait Size % of pages Time: 100 Time: 100 Independent of the pages Multi-page (Landscape) Multi-pages Signal gptions
Destination File setting Printer File prefix: ♥ Screen ♥ Eile ♥ Eile ♥ Unique PDF file name	✓ PDF CSY HIM DIF ✓ IXT XLS EMF
	OK Cancel Help

5 クイックスタートガイド データ分析

Calculate: Percent	Based on:	Area	*		
-ISTD correction					
✓ Use multiplier and dilution fa	actor with ISTDs				
- Calculation factors Use sample data:	From Data File	v #	Compound	ISTD Amount	
Amount	0				
Multiplier:	1				
Dilution:	1				

- 2 上図に示す設定例を使用すると、画面上に面積パーセントレポートを生 成することができます。
- 3 出力先 セクションで、用紙コピー用の プリンタ を選択し、ファイル と PDF を選択して、データファイル内に保存されている PDF レポート ファイルを取得します (.D サフィックスのついたデータファイルは実際 にはディレクトリです。レポートファイルは、ChemStation で直接表示 させるか、あるいは通常の Windows File Explorer を使ってディレクト リ内で見つけることができます)。
- **4** メソッドを再び保存し、レポート設定をメソッドで将来確実に使えるようにします。

メソッドを次に使う場合は、これらの積分イベントとレポートの設定が 使われて、レポートが生成されます。

これで、ChemStation ソフトウェアのデータ分析についての話題は終了します。ChemStation の強力な機能についての詳細は、ChemStation のマニュアルやオンラインヘルプシステムを参照してください。



Agilent 1290 Infinity LC システムマニュアルおよびクイック リファレンス

6 付録

安全に関する情報 114
溶媒情報 117
アジレントのウェブサイト 118
メソッド全体の編集を使ったメソッドの設定 119
メソッド情報 121
機器 / 測定条件 122
データ分析 137
ラン タイム チェックリスト 144

この章では、安全性、法律、ウェブ、およびメソッドの設定に関す る追加情報を記載しています。



安全に関する情報

安全に関する情報

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。

警告と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。

Agilent 1290 Infinity LC システムマニュアルおよびクイックリファレンス

安全に関する情報

ヒューズを交換する際は、必ず指定したタイプ(普通溶断、タイムラグなど)と定格電流のヒューズだけを使用してください。修理したヒューズを 使用したり、ヒューズホルダを短絡させたりしてはなりません。

本書で説明した調整作業には、装置に電源を入れた状態で、保護カバーを 取り外して行うものがあります。その際に、危険な箇所に触れると、感電 事故を起こす可能性があります。

機器に電圧をかけた状態で、カバーを開いて調整、メンテナンス、および 修理を行うことは、できるだけ避けてください。どうしても必要な場合は、 経験のある担当者が感電に十分に注意して実行するようにしてください。 内部サービスまたは調整を行う際は、必ず応急手当てと蘇生術ができる人 を同席させてください。メンテナンスを行うときは、必ず装置の電源を 切って、電源プラグを抜いてください。

本装置は、可燃性ガスや有毒ガスが存在する環境で操作してはなりません。このような環境で電気装置を操作すると、引火や爆発の危険があります。

本装置に代替部品を取り付けたり、本装置を許可なく改造してはなりません。

本装置を電源から切り離しても、装置内のコンデンサはまだ充電されてい る可能性があります。本装置内には、人体に重大な危害を及ぼす高電圧が 存在します。本装置の取り扱い、テスト、および調整の際は十分に注意し てください。

特に、有毒または有害な溶媒を使用する場合は、試薬メーカーによる物質 の取り扱いおよび安全データシートに記載された安全手順(保護眼鏡、安 全手袋、および防護衣の着用など)に従ってください。

安全に関する情報

安全記号

表 10 安全記号

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。
表面が高温の場合に、この記号が装置に付けられます。 加熱されている場合はユーザーはその場所を触れないで ください。

警告は、

警告

注意

人身事故または死に至る状況を警告します。

→ 指示された条件を十分に理解してそれらの条件を満たしてから、 その先に進んでください。

注意

データ損失や機器の損傷を引き起こす状況を警告します。

→ 指示された条件を十分に理解してそれらの条件を満たしてから、その先に進んでください。

溶媒情報

フローセル

ご使用のフローセルの最適な機能を保護するために、以下の基準を守って ください。

• pH 9.5 超のアルカリ性溶液はクォーツに損傷を与え、フローセルの光学 性能を劣化させるため使用を避けてください。

溶媒の使用

溶媒を使用するときは、次の注意に従ってください。

- 褐色の溶媒ボトルを使用すると藻の発生を避けることができます。
- ・ また、次の鉄腐食性溶媒の使用は避けて下さい。
 - ハロゲン化アルカリ化合物およびその酸溶液(ヨウ化リチウム、塩化カリウムなど)。
 - 硫酸や硝酸など高濃度の無機酸(特に高温の場合)。(クロマトグラフィー上可能であれば、ステンレスに対する腐食性の低いリン酸塩またはリン酸緩衝液に変更してください)。
 - ラジカルまたは酸、あるいはその両方を発生するハロゲン化溶媒また は混合液。

 $2CHC1_3 + 0_2 \rightarrow 2COC1_2 + 2HC1$

乾燥クロロホルムを生成する過程で安定化剤のアルコールを除去する と、この反応は速やかに起ります。この反応でステンレスは触媒とし て働きます。

- THF、ジオキサン、ジイソプロピルエーテルなどのクロマトグラフグ レードのエーテルは過酸化物を含む可能性があります。このような エーテルは、過酸化物を吸収する乾性アルミニウム酸化物でろ過して 下さい。
- 強い錯化剤(EDTA など)を含む溶媒。
- 四塩化炭素と 2- プロパノールまたは THF の混合溶液。



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、アジレントのウェブサイトに アクセスしてください。

http://www.agilent.com

Products/Chemical Analysis を選択してください。

このサイトでは、ダウンロード用の Agilent 1200 シリーズモジュールの 最新ファームウェアも提供しています。

メソッド全体の編集を使ったメソッドの設定

1290 Infinity バイナリポンプのすべてのメソッドの手順と設定について 説明します。1290 Infininty クォータナリポンプの場合も、同じ方法で実 施できます。

ChemStation 内のメソッドには、データ取込(システムの制御) 用とデー タ分析(定性的定量的結果を取得するためのデータ処理) 用のすべてのパ ラメータが含まれています。これらのパラメータには、一連の画面を通じ てアクセスできますが、各画面で重点を置いているのは、1 つのモジュー ルまたは機能だけです。これらの画面には、グラフィックユーザーインタ フェース(GUI)のアイコンをクリックするか、メニューバーのドロップダ ウンメニューを使ってアクセスします。新しいメソッドは、既存のメソッ ドまたはブランクのテンプレートメソッド DEF_LC.M を読み込んで編集す ることにより、作成できます。

変更するパラメータの数が少ない場合は、変更対象のパラメータの関連設 定ページに直接アクセスします。慣れていないユーザーの場合は、メソッ ド全体の編集 機能を使う方が簡単です。この機能を使うと、各ページに自 動的にアクセスします。この機能には メソッド > メソッド全体の編集 メ ニューを使ってアクセスします。すると、編集するメソッド項目を選択し てください ダイアログが表示されます。

Edit Method: Instrum	ient 2	
Check Method Section	ns to Edit	
✓ Methy	ed Information	
🗹 įnstru	ment/Acquisition	
✓ Data /	Analysis	
🗹 Bun T	ime Checklist	
ОК	Cancel	Help

図 35 編集するメソッド項目を選択してください

このダイアログには、表示される項目がまとめられており、選択しなかった項目はバイパスできます。

メソッド全体の編集を使ったメソッドの設定

選択した項目に応じて、複数の画面が順番に表示されます。

- ・ メソッド情報は、メソッドに関する説明文で構成されます。
- 機器 / 測定条件 は、以下の項目で構成されます。
 - インジェクタパラメータ
 - ポンプパラメータ
 - オーブンパラメータ
 - 検出器パラメータ
 - 機器カーブ
- ・ データ解析 は、以下の項目で構成されます。
 - シグナル詳細
 - 積分パラメータ
 - レポートパラメータ
- ラン タイム チェックリスト は実行するメソッドの一部を構成します。
- 注記

メソッド全体の編集を実行中に **OK** をクリックすると、現在の入力画面が閉じて次の画面に移動します。これは、一方向プロセスです。

すべてのエントリに入力しないうちに誤って OK を押した場合は、キャンセル を押して メソッド全体の編集 を再起動する必要があります。あるいは作業を 続けていき、最後に未記入の画面に戻るようにします。キャンセル をクリッ クすると、残りの画面を スキップ するためのボタンが表示されます。

メソッド全体の編集を使ったメソッドの設定

メソッド情報

メソッド情報 画面も、メニュー **メソッド** > **メソッド情報** から直接、またはグラフィカルユーザーインタフェースで右クリックすることでアクセスできます。

このボックスには、入力したメソッドについての情報が表示されます。この情報は、メソッドを読み込んでメモリーに常駐させたときに、メソッド **& ランコントロール** 画面のシステムダイアグラムの上に表示されます。

Method Information: Instrument 2	
Method Comments:	
This method runs the test sample to demonstrate how to run an analysis.	<u>s</u>
	×
OK Cancel <u>H</u> elp	

図 36 メソッド情報

メソッド全体の編集を使ったメソッドの設定

機器 / 測定条件

機器メソッドの設定

メソッド セットアップ 画面は、メニュー 機器 > 機器メソッドの設定 から直接、または任意のモジュールアイコン上のグラフィカルユーザーイン タフェースで右クリックし、コンテキストメニュー内の メソッド を選択 することでアクセスできます。メソッド全体の編集 の次の段階は、メソッ ド セットアップ 画面です。この画面には、異なるモジュールまたは機能 のために、6 つのタブが用意されています。

以下のタブがあります。

- 高速オートサンプラ (HiP-ALS)
- ・ HiP-ALS インジェクタプログラム
- バイナリポンプ (BinPump)
- ・ 温度調節機能付きカラムコンパートメント (TCC)
- ダイオードアレイ検出器 (DAD)
- ・ 機器カーブ

タブ間を移動するには、画面上部のタブ名をクリックします。パラメータ を変更すると、適用 をクリックした場合は即座に機器に送信され、すべて のタブに記入して、OK をクリックした場合はすべてのパラメータがモ ジュールに送信されます。そして、画面を閉じて、次の段階に移動します。

アジレントには、機器モジュールに対するコモン RC.Net ドライバという コンセプトがあります。そのため、パラメータの入力タブは、すべてのコ ントロールプログラム (ChemStation、EZChrom、MassHunter など) で似て います。

例題の分離を実行するには、大部分のメソッドと同様に、必ずしもすべて のパラメータを変更する必要はありませんが、完全性を期するために、以 降の項で説明しておきます。

メソッド全体の編集を使ったメソッドの設定

Setup Method		×
👒 HiP-ALS 🔷 HiP-ALS Injector Program 🚆 BinPump 🚀 TCC 😻 DAD	K Instrument Curves	
	HiP-ALS (G4226A)	
Injection Mode	Advanced	-
Injection volume: 1.00 🛟 µL	Auxiliary	
Chandrad intention	Draw speed: 100.0 📫 µL/min	
 Standard injection Injection with peedle wash 	Eject speed: 100.0 ≑ µL/min	
	Draw position: 0.0 🛟 mm	
	Equilibration time: 2.0 📜 sec	
Needle wash	Sample flush out factor: 5.0 📫 times injection volume	
Mode: Flush Port 👻	Vial/Well bottom sensing	Ξ
Time: 3.0 🗘 sec		
Location:	High throughput	
Repeat: 3 🛟 times	Automatic delay volume reduction	
	Enable overlapped injection	
Stoptime Posttime	When Sample Is Flushed Out	
No Limit	After Period Of Time	
	0.00 🗘 min	
		-
	Injection Cleaning	-
Show timetable graph		
	<u>OK</u> <u>Apply</u> <u>C</u> ancel <u>H</u> elp	

オートサンプラタブ (HiP-ALS)

図 37 [メソッド セットアップ] 画面 - [ハイ パフォーマンス サンプ ラ] タブ

- ・ 注入モード
 - 注入量は、注入対象のボリュームを設定します(例:3 µl)。
 - 標準注入は、外部ニードル洗浄は行われていないことを示します。
 - ニードル洗浄+注入は、潜在的キャリーオーバを低減する場合に使います。これは推奨オプションであり、次のエントリで構成します。
- ・ ニードル洗浄 は、上記で選択した場合です。
 - モードは、ニードル外部を洗い流す方法を指定します。フラッシュ ポートで強力に洗い流すか、指定した洗浄バイアルに浸します。

メソッド全体の編集を使ったメソッドの設定

- フラッシュポートに接続されているペリスタルポンプが、洗浄溶媒を
 送り出す 時間 を秒単位で指定します。ペリスタルポンプは、フラッシュポートを洗浄するために、さらに 15 s 送り出します。
- ロケーションは、洗浄バイアルを選択している場合に、バイアルまたはウェルプレートのいずれを使うのかを指定します。
- 注記 セプタムを持つバイアルは使えません。すなわち、セプタム上のキャリーオー バ剤が輸送されることを回避するために、オープンである必要があります。
 - 繰り返しは、洗浄バイアルを選択している場合に、ニードルをバイアルに浸す回数を指定します(デフォルトは3、最大で5)。
 - ストップタイム/ポストタイム には 無制限/オフ が設定されますが、ポンプタブではこれらの値は慎重に取り扱う必要があります。
 - 詳細設定 補助設定
 - 吸引速度は、サンプルがニードルに吸い込まれる速度です。デフォルト値は、100 µl/minです。粘性のあるサンプルの場合は、少ないサンプル量(<2 µL)でも十分な精度を確保するために速度を落とす必要があります。
 - 吐出速度は、ニードルから吐き出す速度です。
 - 吸引ポジションは、バイアル底部の上の標準注入ポジション10 mm からの垂直オフセットです。これは2mlバイアルの約半分なので、 サンプルをバイアル底部近くから取り出すには負のオフセットが必要 になります。たとえば、値-7 mmを指定すれば、ニードル3 mmの チップがバイアル底部の上に配置されます。
 - 平衡化時間は、サンプルを吸い込んでからニードルを移動させるま
 での間のディレイタイムです。
 - サンプルフラッシュアウト係数は、注入後にバルブがバイパスに切り替わるまでオートサンプラが待機する時間を決定します。この係数によって、ニードル、シート、注入バルブを使ってサンプルゾーンから汚れが取り除かれることが保証されます。デフォルト値は5です。
 - バイアル/ウェル底部センサは、吸引ポジションのオフセットの使用の代替です。ニードルは、バイアルまたはウェルの底部に接触するまでゆっくりと下降し、その後1mmだけ上昇します。これは、ニードルをバイアル底部に近付けることを保証するための優れた方法ですが、注入の完了までに時間がかかることと、バイアル底部に粒子状物

メソッド全体の編集を使ったメソッドの設定

質がある場合はニードルを詰まらせる可能性があるため、ニードルが 使えなくなるという欠点を持ちます。

- ・ 詳細設定 ハイスループット
 - 自動バルブ切り替え(ADVR)は、注入が完了し、サンプルフラッシュ アウト係数で定義されたボリュームがインジェクタを通過した後に、 注入バルブをメインパスからバイパスへ切り換えます。この係数に よって、システムディレイボリュームは70 µ1 ほど削減され、グラ ジエント変更がカラムに到達する時間が短縮されます。
 - オーバーラップ インジェクション も、サンプルがインジェクタから フラッシュアウトされた後か、または分析中に指定されてている時間 が経過した後のいずれかの時点で、注入が完了すれば、注入バルブを メインパスからバイパスへ切り換えます。インジェクタは次回の注入 に備えるために次のサンプルを吸引するので、サイクルタイム全体を 短縮し、サンプル処理数を増加させます。
- 注入バルブクリーニング
 - インジェクタクリーニングを使うと、注入システムを溶媒でフラッシュできます。
 - 注入バルブクリーニングを使うと、対象化合物が注入されたときに キャリーオーバを最小にするため、分析中の設定値でバルブを切り換 えられます。

メソッド全体の編集を使ったメソッドの設定

ハイ パフォーマンス サンプラ (Hip_ALS インジェクタ プログラ ム) タブ

Setup Method	X
👒 HiP-ALS 🧇 HiP-ALS Injector Program 🚆 BinPump 🚀	TCC 🔝 DAD 🦗 Instrument Curves
Use Injector Program	
Function	Parameter
Draw -	Draw default volume from sample with default speed using default offset
Append Insert Delete Clear all	Move <u>up</u>
Cut Copy Paste	Move <u>d</u> own
Show timetable graph	
	<u>UK</u> <u>Apply</u> <u>Cancel Help</u>

図 38 [メソッド セットアップ] 画面 - [HiP サンプラ インジェクタ プ ログラム] タブ

これを使うと、たとえば、プレカラム誘導体のような複数のバイアルのア リコートの操作を含む、特殊な注入手順を作成できます。検出機能または 感度を向上させるために、サンプルには試薬が自動的に混合されます。よ く使われる例は、OPA および FMOC 試薬を用いるアミノ酸誘導体です。詳 細は、Agilent 1290 Infinity Autosampler manual を参照してください。

メソッド全体の編集を使ったメソッドの設定

ietup Method							X
🧇 HiP-ALS 🧇 HiP-ALS Injector Program	🚆 BinPump 🚀 TCC 🔻 DAD	🔀 Instrument Curve	s				
					BinPum	p (G4220A)	
Flow		Timetable					A
0.	000 🛟 mL/min	Time 🗠	Function			Parameter	
Solvents							
1 @ H20	•						
2 © H20	•						
B: 00 1 %	•						
2 O H20	•						
Stoptime	Posttime						
No Limit	Off						
1.00 ; min	1.00 ; min						
Property Limite							
Nin 0.00 the	Mars (00.00 t) has	Add	<u>R</u> emove	Clear all			
Miri. U.UU Jar	Max. 400.00 Dar	Cut	Сору	Paste			
							-
Show timetable graph							

バイナリポンプ (BinPump) タブ

図 39 [メソッド セットアップ] 画面 - [バイナリ ポンプ] タブ

- 流量には、最大で 5 ml/min の流量が設定されます。たとえば、分離
 0.4 ml/min が使われます。背圧が最大圧力設定値に簡単に到達すれば、流量は数秒間減少し圧力を低下させますが、圧力がこの方法で制限され続けると、エラー状態が生成されて流れが停止します。
- 溶媒 は、利用可能な移動相と 2 つのチャンネル A と B に送り出す割 合を定義します。各チャンネルのドロップダウンボックスを使えばリス トから溶媒を選択できるので、ポンプコントロールで最適の圧縮率設定 を使えるようになります。これにより、『「最適ディレイボリュームの設 定方法」42ページ 図』で説明したように、フロー特性を最適化できま す。2 番目のテキストボックスには、入力する移動相についての説明を 記述します。ポンプに溶媒切り替えバルブが取り付けられている場合は 各チャンネルは 2 つの溶媒オプションを持つことになり、溶媒の説明文 の左にあるラジオボタンを使えばメソッドに適したオプションを選択で きます。ポンプは指定された A および B チャンネル (たとえば、A2 と

メソッド全体の編集を使ったメソッドの設定

B1) のバイナリ混合を行います。A1 と A2 または B1 と B2 の混合は許 されていません。A と B の割合として入力される値は、アイソクラ ティックメソッドの混合、またはグラジエントメソッドの開始条件とグ ラジエント分析間の平衡化条件を定義します。値は B についてのみ入力 します。カーソルを移動させれば、A には 100% - B の値が自動的に入 力されます。例題の分離に対して、A には水、B にはメタノールを 50 % の割合で設定します。A は 50 % になります。

• タイムテーブル には、移動相内の A と B の混合の割合、または、必要 に応じて、許される流量と最大圧力を使って、分析中に生ずる変化の詳 細が設定されます。タイムテーブルは、指定した設定値間でパラメータ をリニアに変化させます。この画面の別の場所で指定する設定は初期状 態として使われ、変更されるのは、タイムテーブルにエントリを作成し たときだけです。たとえば、分析中は流量が一定ならば、タイムテーブ ルに流量に関するエントリを作成する必要はありません。タイムテーブ ルにエントリを作成するには、以下の手順を実行します。追加 ボタン をクリックして、タイムテーブルに行を追加します。設定値の時間を入 力します。ドロップダウンリスト(混合、流量圧力)からエントリのタ イプを選択します。そして **パラメータ** ボックスをクリックして、値の エントリボックスを表示します。タイムテーブルのエントリを論理シー ケンスで作成した場合は、エントリは時刻に従って、自動的に順番が並 べ替えられます。タイムテーブルの行は直接編集できます。また行の追 加と削除には、切り取り、貼り付け、削除 などのボタンを使用できま す。グラジエントプロファイルを作成するために、一連のリニアグラジ エントセグメントを指定するための複数の行を追加できます。例題の分 離に対して単純なグラジエントを設定するには、まず、すでに空になっ ていなければ、全消去 ボタンを使ってタイムテーブルをクリアし、 4.00 min に行を追加し、混合を 90 % に変更します。 グラフ画面には、 50 % B から 90 % B への 4 min 間にまたがるリニアグラジエントが表 示されます。ステップグラジエントが必要な場合は、2 つのエントリ 「before step」と「after step」の設定(0.01 min で区切る)を追加す ることで作成できます。これは、分析の最終段階で、ステップ グラジエ ント内で強力な溶媒や流量を増やす(たとえば、%Bを75%から 95%にステップアップする)ことで、強力に保留されているピークを 短時間でカラムから溶出させる場合に使われます。タイムテーブル内に 時間設定 0.00 min を入力する必要はありません。この値は、この画面 の別の設定値から持ってこられます。ただし、タイムテーブル内の完全 なリストを表示させる場合は、0.00 min のエントリを作成します。これ については何の問題もありませんが、初期状態が変更された場合は、新

128

メソッド全体の編集を使ったメソッドの設定

しい設定をタイムテーブルと、画面の溶媒セクションの設定値の両方に 入力する必要が生じます。

- タイムテーブルグラフの表示は、ボックスにチェックが付けられた場合は、タイムテーブルの変化をグラフで表示します。
- ストップタイム は、分離または分析の全体の時間を定義します。「ラン タイム」と呼ばれることもあります。これは注入が行われてから分析が 終了する(つまり、データ取込は停止し、フロー、混合、その他のシス テム設定はメソッドの初期値に戻され、システムは次の注入に備えて準 備が終わる)までを分単位でカウントする時間です。この値は、タイム テーブルの最後のエントリの値以上である必要があります。そうしない と、タイムテーブルイベントが完了する前に、分析が停止して開始条件 に戻ってしまいます。ストップタイム には 無制限 を設定できます。こ の場合は、分析を手操作で停止する必要があります。すべてのシステム モジュールはストップタイムパラメータを持っていますが、ポンプのス トップタイムがマスタと見なされます。そのため、他のモジュールは、 通常、この値に従って設定されます。
- ポストタイム は分析終了後のカウントダウン時間を定義します。この時間内は次の注入が禁止されます。これは、システムがグラジエント分析後に再平衡化するための時間です。アイソクラティックメソッドの場合は、オフ を設定します。グラジエントメソッドの場合は、値はベースライン処理を観察することによって経験的に決定されますが、通常はシステムのディレイボリュームに、システムでフラッシュされるカラムボリュームを少なくとも 3~5 つ加えた時間が必要になります。

メソッド全体の編集を使ったメソッドの設定

Setup Method			×
👒 HiP-ALS 👒 HiP-ALS Injector Program 🕌	BinPump 🚀 TCC 🔝 DAD	M Instrument Curves	
			TCC (G1316C)
Temperature		Timetable	
Left	Right:	Advanced	<u>^</u>
Not Controlled	Not Controlled	Enable Analysis	
● <u>40.0</u> *C	D 20.0 📜 *C	when front door open	Dista
As Detector Cell	As Detector Cell	Lerc	night.
	Combined	With any temperature	With any temperature
Column Switching Valve		When temperature is within + 0.8 * *C	When temperature is within + 0.8 ⁺ °C
Port 1 → 2			
Stoptime Postti	me		
As Pump/Injector As Pump/Injector I.00 i min	● Off ● 1.00 ; min		*
Show timetable graph			
			ply <u>C</u> ancel <u>H</u> elp

温度調節機能付きカラムコンパートメント (TCC) タブ

図 40 [メソッド セットアップ] 画面 - [カラムコンパートメント] タ ブ

 ・ 温度 は、左側と右側のカラムホルダの温度を定義します。独立に制御 することも、結合 ラジオボタンをオンにして同時に制御することも可 能です。結合した場合は、左側の設定によって両方のセクションが制御 されます。これが必ず必要になるのは、カラムが 15 cm よりも長く、両 方のセクションでサポートが必要な場合です。2 つのカラムが異なる温 度で動作する必要がある場合は、両側は独立に運用します。このような 状況は、カラム切り換えバルブが実装されている場合に発生します。独 立した温度ゾーンを使う別の例としては、カラムが一方のサイドでは高 温(たとえば、60 ℃以上)で動作しており、他のサイドの熱交換器が、 検出器に渡す前にフローセル内の温度効果が原因のノイズを減らすため に溶離液の冷却で使われる場合です。検出器セルと同じ オプションを選 択すると、自動的に、1290 Infinity 検出器内のセルから温度が読み取 られます。

メソッド全体の編集を使ったメソッドの設定

各ゾーンの温度には、-5 ℃ ~ 100 ℃ の範囲で設定できますが、カラ ムがその温度で動作可能なことを確認しておく必要があります。(範囲 の上端では、Agilent ZORBAX RRHD および RRHT StableBond 相が使えま す。)温度は周囲温度の ± 0.15 ℃ ~ 10 ℃ の範囲になるように制御 されますが、12 ~ 15 ℃以下の温度で動作するアプリケーションはほと んどないことに注意してください。このような低温で TCC を使うことは 避けてください。湿気を含んだ空気が凝縮して水になり、リークセンサ をトリガするからです。

- カラムスイッチングバルブは、バルブがカラムホルダに適合する場合 だけアクティブになるオプションです。次の3種類のバルブが用意され ています。
 - 2 ポジション 6 ポート 2 カラム間を切り換えるために使用
 - ・ 2 ポジション 10 ポート カラム再生を切り換えるために使用
 - 8 ポジション 9 ポート MDS 内で複数のカラムを選択するために使用

メソッド全体の編集を使ったメソッドの設定

Setup Method	🔀
👒 HiP-ALS 👒 HiP-ALS Injector Program 🞽 BinPump 🖋 TCC 🖤 DAD	X Instrument Curves
	DAD (G4212A)
Signals	Timetable
	Advanced
Use Wave Band Reference Reference Signal length width Wavelength Bandwidth	Spectrum
	Store: All
Signal A 🛛 250.0 ; 100.0 ; 🖓 360.0 ; 100.0 ; nm	
Signal B 🖸 254.0 ; 16.0 ; 🖸 360.0 ; 100.0 ; nm	Range from: 190.0 🛟 to 400.0 🛟 nm
Signal C 210.0 7 8.0 7 360.0 7 100.0 7 nm	Step: 2.0 ⁺ nm
Signal D 230.0 16.0 1360.0 100.0 mm	
Signal E 280.0 16.0 1360.0 100.0 mm	Threshold: 10.0 , mAU
	Analog Output
Signal H = 2000 1 100 1 = 3600 1 1000 1 mm	Output 1:
	Zero Diffeet: 5 * %
Peakwidth	
>0.025 min (0.5 s response time) (10 Hz)	Attenuation: 1000 • mAU
	Margin for negative Absorbance Slit
Stoptime Posttime	
As Pump/Injector Off	Autobalance Lamps on required for analysis
	V Prerun V UV Lamp
Show timetable graph	
	<u>O</u> K <u>Apply</u> <u>Cancel Help</u>

ダイオードアレイ検出器(DAD)タブ

図 41 [メソッド セットアップ]画面 - [ダイオードアレイ検出器]タ ブ

- ・ シグナル:最大で 8 つの独立したシグナル (クロマトグラム)を記録で きます。シグナルに収集とマークするには、そのシグナルに対応した 使 用するシグナル ボックスにチェックを付け、波長とバンド幅を定義し、 リファレンスシグナルが必要な場合は、ボックスにチェックを付けて定 義します。
 - 波長には、シグナルの中心波長(nm)を設定します。
 - バンド幅 には、シグナルの幅 (nm) を設定します。
 - リファレンス波長 には、分析シグナルから抽出したリファレンスバンドの中心波長 (nm)を設定します。
 - リファレンスバンド幅には、リファレンスバンドの幅(nm)を設定します。

メソッド全体の編集を使ったメソッドの設定

- ピーク幅には、データ収集レートとシグナルフィルタリングを設定します。
- ストップタイム / ポストタイム には 無制限 / オフ を設定します。こ れらの値は、通常、ポンプ タブで取り扱います。ポンプで定義された分 析の終了よりも前にデータ分析を停止する必要がある場合は、検出器モ ジュールのストップタイムが、ポンプのストップタイムと異なることが あります。グラジエントの終了時にグラジエント平衡化ランプが設定さ れる場合がこれに相当します。たとえば、%B が 95 % 95 % にまで上昇 (10分分後)し、その間に、すべてのピークがカラムから溶出すると します。この場合、分析は本質的には終了していますが、エクストラグ ラジエントセグメントが追加されているため、2 分間にわたって初期値 は %B 後退します。この間にカラムの再平衡化がなだらかに開始されま す。この下方ランプでは有益なデータは期待できないので、検出器のス トップタイムには 10 分 を設定しデータ収集を停止します。一方、ポン プのストップタイムには 12 分 を設定して下方ランプが完了するのを待 ちます。またユーザーによっては、検出器にだけ異なるストップタイム を設定するのが煩わしいため、クロマトグラムの最後の数分には有用な データを含められないことを納得した上で、そのまま記録させることを 選択する場合もあります。検出器のストップタイムは、他のモジュール の早期のストップタイムと同様に、ポンプのストップタイムには影響を 与えずに、分析を終了させます。そのため、ストップタイムに ポンプ 同様 という設定が用意されています。
- タイムテーブルは、他のモジュールと同様に操作します。すなわち、 行を追加し、変更する機能を選択し、その機能に新しい値を設定します。検出器の変更は、指定に伴って直ちに行われます。以下の機能は、 分析中に変更できます。
 - バランス
 - シグナル変更
 - しきい値変更
 - ピーク変更 検出器のピーク幅
 - スペクトル取得モードの変更
 - 接点の変更

・ 詳細設定 - スペクトル

スペクトルは、分析中に、連続的またはピーク制御ベースで保存されま す(これは ChemStation ソフトウェアの場合です。たとえば、EZChrom

133

メソッド全体の編集を使ったメソッドの設定

のようなソフトウェアは、すべてのスペクトルについて連続的な収集し かサポートしていません。ピーク制御オプションは表示されません)。 スペクトル収集とシグナル収集は、検出器のファームウェアが実行する 独立した操作であり、コンピュータのソフトウェアが 3D データマト リックスから抽出するデータとは無関係です。スペクトルを収集する レートは、ピーク幅の設定で指定します。ピーク幅で指定した時間内 に 8 つのスペクトルが収集されます。ピーク制御スペクトルを保存する 場合は、ファームウェアが実行するピーク検出はシグナル A だけが対象 です。シグナルが複数ある場合は、シグナル A をブロードバンド検出器 として設定し、波長の異なるすべてのシグナルについてピークスペクト ルを収集できるようにする必要があります。

保存は、以下のオプションを使って、スペクトル収集モードを制御します。

なし - 保存するスペクトルはありません。

頂点 + ベースライン - ピークの開始点、頂点、終了点の 3 つのスペ クトルを収集します。

頂点+スロープ+ベースライン - ピークの開始点、アップスロープ、 頂点、ダウンスロープ、終了点の 5 つのスペクトルを収集します。

ピーク内全て - ピーク内の全てのスペクトルが保存されます。

全て - 分析を通じて全てのスペクトルが保存されます。

2 スペクトル毎 - 分析を通じて取得したスペクトルが 1 つおきに保存されます。

- 範囲 検出器の 190 nm ~ 640 nm の全範囲にわたって、またはユー ザーが指定した限定された範囲にわたって、スペクトルが保存されま す。これにより、保存するデータ数を減らすことができます。
- ステップ スペクトルに保存するデータの間隔 (nm)、つまりスペクトルの解像度を制御します。大部分のアプリケーションにおいて、デフォルト設定 2 nm が優れた選択です。
- スレッショルド この値(mAU)よりも低いピークしか持たないスペクトルは保存しません。

詳細設定 - アナログ出力

1290 Infinity 検出器は、デジタル出力を受け付けないデータシステム 用にアナログシグナル出力コネクタを 1 つ持っています。以下のパラ メータを設定できます。

メソッド全体の編集を使ったメソッドの設定

- ゼロオフセットは、出力シグナルに設定する割合として、ゼロレベルを設定します。これにより、負のドリフトを持つスコープでも処理できるようになります。
- アッテネーションは、設定された吸収率がフル出力となるようにスケーリングします。
- ・ 詳細設定 ネガティブ吸収のマージン

デフォルト設定は、100 mAUです。これは、検出器が十分なダイナミックレンジを持つことを意味します。ゼロレベルが設定されている場合でも、この値に合わせて計測できます。大きな負のピークを計測するには、または大きな負のドリフトを持つベースラインに従うには、レンジの底部でシグナルが平らになってしまうことを避けるために、この値を下方修正する必要があります。ただし、正当な理由がない限り、変更してはいけません。大きな負の値を受け入れるようにすると、ベースラインノイズが増加し、正のピークを計測するレンジが減少してしまうからです。

• 詳細設定 - スリット

分光器への入射スリットは、スペクトルの解像度を制御し、ベースラインのノイズと感度に影響します。デフォルト設定は、4 nm ですが、これは大部分のアプリケーションに適しています。このパラメータについての詳細は、『「感度を向上させる方法」59 ページ 図』を参照してください。

- 詳細設定 自動バランス は、すべての波長に対して吸光レベルにゼロ を設定します(つまり、スペクトルのすべてのポイントをゼロと平衡化 させます)。これにより、ベースラインシグナルもゼロになります。プ レラン は、分析を開始する直前に平衡化させるために選択します。こ れが一般的な選択です。代わりに、ポストランタイムが経過した後、分 析の終了時に平衡化するために、ポストラン が選択されることもあり ます。たとえば、シグナルが常に負のドリフトを示しているため、分析 をゼロの吸光率で終了させたいと考えた場合は、このパラメータを設定 することによって、次回の分析で正しいセロレベルが設定できます。こ のパラメータは、分析の終了時に行った平衡化を、過去に遡って変更す るわけではありません。
- ・ 詳細設定 分析時のみランプをオンにする: 1290 Infinity DAD または
 MWD は、重水素ランプを 1 つ持っています。分析で必要になる場合は、
 このボックスにチェックを付ける必要があります。

メソッド全体の編集を使ったメソッドの設定

機器カーブ タブ

ietup Method
🧇 HIP-ALS 🚳 HIP-ALS Injector Program 🕌 BinPump 🚀 TCC 🔝 DAD 🧮 Instrument Curves
BinPump:
✓ Pressure (bar)
✓ Flow (ml/min)
✓ Solvent Ratio A (%)
✓ Solvent Ratio B (%)
TCC:
Left Temperature (*C)
Right Temperature (*C)
DAD:
Board Temperature (*C)
Dptical Unit Temperature ("C)
UV Lamp Anode Voltage (V)
Show timetable graph
<u>OK</u> <u>Apply</u> <u>Cancel Help</u>

図 42 [メソッド セットアップ] 画面 - [機器カーブ] タブ

[機器カーブ]タブを使うと、関連ボックスにチェックを付けることに よって、検出器シグナル以外の監視対象データストリームをデータと共に 保存できます。これは、主として診断目的で使われます。以下のタブがあ ります。

- ・ ポンプ:
 - 圧力
 - 流量
 - A/B 合成 クロマトグラムにグラジエントプロファイルを重ね合わせるために役立ちます。
- 温度調節機能付きカラムコンパートメント
 - 左側 / 右側温度
- 検出器:
 - ボード温度
 - オプティカルユニット温度
 - UV ランプ陰極電圧

メソッド全体の編集を使ったメソッドの設定

データ分析

シグナル詳細

メソッド & ランコントロール ビューからは、シグナル詳細 画面にも直接 アクセスできます。キャリブレーション アイコン上の GUI を右クリック して、コンテキストメニュー内の シグナル詳細 を選択します。[データ分 析] ビューでは、キャリブレーション > シグナル詳細 を使ってアクセス できます。

シグナル詳細 画面は、メソッド全体の編集 の次のステップであり、取得 したシグナルのうち、処理対象とするシグナルを [データ分析] に対して 通知します。ドロップダウンボックスには、検出器設定で定義された分析 シグナルと、温度、流量、混合、圧力、診断トレースのような記録された パラメータを含む、利用可能なシグナルがリストされます。シグナルを選 択し メソッドに追加 をクリックして、シグナルを画面下部にある シグナ ル詳細 テーブルに送信します。取得した検出器シグナルの一部またはすべ てを処理対象として選択できます。シグナルを一切選択しなかった場合は テーブルは空です。この場合は、ChemStation はデフォルトとして、取得 した検出器シグナルのすべてを処理対象とします。

既存のメソッドを編集して新しいメソッドを作成した場合は、メソッドの 実行を試みると、パラメータ ミスマッチ エラーが表示されることがあり ます。この理由は、旧メソッドでは シグナル詳細 に、特定のシグナル、 たとえば、250 nm 、8 nm バンド幅が含まれていたけれども、新メソッド では、それを、たとえば、254 nm/12 nm に変更してしまったからです。シ グナル詳細 テーブルには、元のシグナルが残っており、取得されることが なくなったシグナルを処理するというメッセージが表示されます。テーブ ル内の元のシグナルを強調表示し、行削除 ボタンを押して問題を解決しま す。

システムが、ダイオードアレイ検出器、質量分析計のように、複数の検出 器を使っている場合は、シグナル説明の行を使って、下流の検出器のディ レイ時間を入力し、ソフトウェアが異なる検出器からのピークを調整でき るようにします。

137

メソッド全体の編集を使ったメソッドの設定

Signal Details: Instrument 2		$\overline{\mathbf{X}}$
Available Signals DAD1 A, Sig=250,100 Ref=off	Add to Method	
Insert Row Append Row Delete Row Signal Description	Start End Delay Alig	ın Peak 1 Peak 2 Align Window
	OK Cancel <u>H</u> elp	

図 43 シグナル詳細

メソッド全体の編集を使ったメソッドの設定

積分イベントの変更

[積分イベント] アイコン上のグラフィック ユーザー インターフェース を右クリックして、コンテキストメニューの 積分イベントの編集 をク リックすることにより、メソッド & ランコントロール ビューからは、積 分イベントの編集 画面にも直接アクセスできます。データ分析 ビューで は、積分 > 積分イベント... メニュー、または 積分イベントの編集 タス クアイコンを使ってアクセスできます。

Single Runs: PHENONES	
<u> Integration</u> 🎂 Calibration 📶 Sig	gnal 🛄 P
📠 1 Keport: Short	1
DAD1 A, Sig=242ONE-MIX-008.D) 🔻 🎳	Ð, Ə,
> ₹ 3 0 0	

Method Manual Events

Initial Events For All Signals:

Integration Events	Value
Tangent Skim Mode	Standard
Tail Peak Skim Height Ratio	0.00
Front Peak Skim Height Ratio	0.00
Skim Valley Ratio	20.00
Baseline Correction	Classical
Peak to Valley Ratio	500.00

Specific Events For Signal:

DAD Default 👻

Tim	e	Integration Events	Value
Init	al	Slope Sensitivity	5
Init	al	Peak Width	0.05
Init	al	Area Reject	5
Init	al	Height Reject	1
Init	al	Shoulders	OFF

図 44 [積分イベントの編集] 画面

積分、キャリブレーション、レポーティングは、メソッドのデータ分析部 分です。積分パラメータとキャリブレーションテーブルは、データを取得 し、データを[データ分析]ビューでレビューすることで簡単に設定でき ます。積分イベントはその時点で最適化できます。また、最初の取得の分 析では、通常、デフォルトの設定を使います。

積分イベントの編集 画面には、2 つのテーブルがあります。

 全てのシグナルの初期イベントは、メソッドで取得するすべてのシグ ナルに適用されるイベント(積分パラメータ)を含みます。

139

メソッド全体の編集を使ったメソッドの設定

 ・ シグナルの特定のイベントは、1 種類の検出器または同一検出器からの 異なるシグナルに固有のイベントを含みます。

このテーブル内の重要なパラメータは、以下のとおりです。

- スロープ感度は、ピークの開始と終了をマークするのに必要な、
 ベースラインのスロープと曲率を表わします。
- ピーク幅:対象のピークの中で最も狭いピークの半分の高さの幅を設定します。これは、インテグレータが、ノイズと非常に小さなピークを区別することを助けます。
- 面積リジェクト / 高さリジェクト は、棄却を制御する値です。この値より小さい面積または高さは、ピークの結果からは棄却されます。
- 積分オフ/オンは、設定されたリミット間の積分を抑止します。ほとんどの場合、注入から、溶媒の前または保留されていないピークマーカーまでの範囲の積分を禁止するために使われます。

積分オフ / オンのような行は、ウィンドウの最上部にあるアイコンを使って、テーブルに追加します。

OK をクリックして終了すると、[メソッド全体の編集] プロセス内に次の 画面が表示されます。

メソッド全体の編集を使ったメソッドの設定

レポート条件

[レポート] アイコン上のグラフィック ユーザー インターフェースを右 クリックして、コンテキストメニューの レポート条件 を選択することに より、メソッド & ランコントロール ビューからは、レポート条件 画面に も直接アクセスできます。データ分析 ビューでは、レポート > レポート 条件 メニュー、または [レポート条件] タスクアイコンを使ってアクセ スできます。

Specify Report: Instrument 1				×
Reporting settings Quantitation settings				
Report mode ○ Use Intelligent Reporting	c Reporting			
Style Report style: Short Quantitative result sorted by: Signal	× ×		 Chromatogram output Pogtrait 	Size % of pages
 ☐ Sample info on each page ✓ ≜dd chromatogram output ☐ Add sample custom fields to sample info 	✓ Add fraction table and ticks Add summed peaks table Add compound custom fields		 Landscape Multi-page (Landscape) 	Response: 40
Report layout for uncalibrated peaks O Separately O With calibrated peaks	O Do not report		1 🗘 Pages	Signal <u>o</u> ptions
Destination File setting □ Drinter File prefix Report ☑ Screen Prefix Report	♥ PDF	□ CS <u>V</u>	<u>н</u> тм	DIF
V Eile	<u>⊾</u> <u>T</u> ×1	<u>∏ ×</u> LS	<u> </u>	
			ОК	Cancel Help

図 45 [レポート条件] 画面

プリンタや PDF ファイルに出力するためのクラシックレポートによる単純 な 面積% レポートを設定するには、[レポート条件] 画面の以下のセク ションに以下の設定を入力します。

レポート設定 タブ:

- レポートモード: クラシックレポートを使用する
- ・スタイル
 - レポート スタイル: 簡易
 - ・ 定量結果の表示順: シグナル

メソッド全体の編集を使ったメソッドの設定

- クロマトグラム出力の追加: チェックあり
- クロマトグラム出力:縦
- サイズ:
 - 時間軸 100 %/ ページ
 - レスポンス軸 40 %/ページ
- ・ 出力先
 - プリンタ: チェックあり
 - 画面: チェックなし
 - **ファイル**: チェックあり
- ファイル設定:
 - **PDF**: チェックあり
 - 固有 PDF ファイル名: チェックあり

定量設定タブ:

- ・ 計算モード
 - 計算:パーセント
 - カウント法: 面積

OK をクリックして終了すると、[メソッド全体の編集] プロセス内に次の 画面が表示されます。

メソッド全体の編集を使ったメソッドの設定

機器カーブ

Instru	nent	Curves: Instrument 2	×
C 1			
Selec	t curv	es to overlay:	
		Instrument Data Curves:	
		%A 🔺	
		%B	
		%C(C1)	
		%D(C2)	
		Flow	
		Temperature	
		High Pressure	
		MSD: Ion Polarity	
		MSD: Polarity - 1	
0	К	Cancel <u>H</u> elp	

図 46 [機器カーブ] 画面

機器カーブ チェックボックスを使うと、クロマトグラムに記録されたパラ メータをグラフとして重ねることができます。 メソッド全体の編集を使ったメソッドの設定

ラン タイム チェックリスト

ラン タイム チェックリスト 画面も、メソッド > ラン タイム チェック リスト メニューから直接、または画面の右上の ラン タイム チェックリ スト アイコンをクリックすることでアクセスできます。

Run Time Checklist: Instrument 2		×
Check Method Sections to Run		
Pre-Run Command / Macro		
✓ Data <u>A</u> cquisition		
☑ Standard <u>D</u> ata Analysis		
Customized Data Analysis Macro		
Save GLP Data		
Post- <u>B</u> un Command / Macro		
Save <u>M</u> ethod with Data		
OK	Cancel Help	

図 47 [ラン タイム チェックリスト]画面

ラン タイム チェックリスト は、メソッドでデータ収集とデータ分析の両 方を実行して、マクロコマンドまたはプログラムをワークフローの各種の ポイントにリンクさせるか否かを選択します。多くの場合、データ収集 と 標準データ解析 チェックボックスにチェックを付けます。たとえば、一連 のメソッドを開発しているときのように、データ分析が不要の場合は、標 準データ解析 はチェックを外して、レポートを生成しないようにできま す。データは後から データ分析 ビューを使って表示できます。

マクロプログラム内でアクセスポイントの 1 つにリンクするには、関連 ボックスにチェックを付け、右にあるテキストボックスにマクロ名を入力 します。ソフトウェアは、C:¥Chem32¥Core ディレクトリ内のマクロを検索 します。他の場所にある場合は、パスを含めます。

メソッドのワークフロー内のアクセスポイントは、以下のとおりです。

・ プレランコマンド / マクロ
付録 6

メソッド全体の編集を使ったメソッドの設定

- カスタマイズされたデータ分析マクロ
- ポストランコマンド / マクロ

データファイルにメソッドを保存は、メソッドのコピーをデータファイル RUN.M に保存します。ChemStation を通常の構成で実行している場合は、 この作業は不要です。このソフトウェアによって、このデータファイルが 常に保存されているからです(B.02.01 以降のすべてのバージョン)。この オプションが意味を持つのは、たとえば、ChemStation で、**固有シーケン** スフォルダの作成 をオフにし、メソッドが自動的にデータファイルにコ ピーされないようにする設定を行っている場合だけです。

以上のプロセスの中でこの画面が最後の画面です。OK をクリックすると、 ラン タイム チェックリスト が終了し、メソッド全体の編集 プロセスが 終了します。この段階で、ファイル > 名前を付けて保存 > メソッド また は メソッドメニュー > 名前を付けてメソッド保存を使うと、メソッドは マスターメソッドディレクトリ (デフォルトでは、C:¥Chem32¥1¥Methods) に保存されます。

索引

A

Agilent 1290 Infinity LC シ ステム システムコンポーネン ト 25 新機能 22 性能の範囲 22

D

DEF_LC.M 119 I Infinity オートサンプ ラ 説明 32

J

Jet Weaver 説明 48

۷

van Deemter 方程式

ア

アジレント インターネット上 118 安全 一般的な情報 114 安全クラス I 114 安全 記号 116 <mark>イ</mark> インターネット 118

え

液体クロマトグラフィ より小さな粒子を使用す る 8

か

カラム コンパートメン ト 説明 34 カラム サブ 2 マイクロンの粒 子 14 カラム外ボリューム 40 カルキュレータ コスト 14

き

機器 / 測定条件 120

<

9

クイックスタートガイド はじめに 92

け

検出器 高感度の達成 61

E

コンフィグレーション

1 スタック 74, 79 2 スタック 77, 82 2 スタック前面 77, 82 2 スタック背面 78, 83

さ

最適化
HPLC の条件 11
クロマトグラフィの分離性
能 11
検出器感度 61
スリット幅 64
注入量 51
波長と帯域幅 61
サブ 2 マイクロンの粒
子 14

L

シグナル 積分 109 シグナル波長 63 システムコンポーネント オートサンプラ 32 カラム コンパートメン ト 34 ダイオードアレイ検出 器 36 バイナリポンプ 25

す

スリット幅 64

146

せ

積分 シグナル 109

た

ダイオードアレイ検出器 説明 36

ち

注入量		
ボリ	ュームを増加させ	
る	51	

で

ディレイボリューム		40
データ解析	120	
データ取込レ-	- ト	65

は

バイナリポンプ 説明 25 バンド幅 63 波長と帯域幅 最適化 61

v

ピーク幅 65

ふ

フローセル Max-Light カートリッジフ ローセル 61 Max-Light 高感度セ ル 61 分解能 11 分離能 17

ほ

保持係数 11 ポンプのパージ 87

め

```
メソッド情報 120
メソッド
デフォルト 119
メソッド全体の編
集 119
```

6

ランタイムチェックリス ト 120

れ

レスポンスタイム 65 レポート 仕様 111

ろ

理論段数 11

www.agilent.com

本書の内容

本書には、Agilent 1290 Infinity LC システ ムの技術的リファレンス情報が記載されてい ます。

本書では、以下の項目について説明します。

- 概要
- 製品の説明
- ・ システムの最適化
- 設定と設置
- クイックスタートガイド

© Agilent Technologies 2009-2011, 2012

Printed in Germany 05/2012



G4220-96301

