医薬品



# 糖ペプチドベースのキラルカラムによる 最新のキラル分離

Agilent InfinityLab Poroshell 120 キラル-T カラムを用いた マレイン酸チモロールのキラル分離

#### 著者

William J. Long Agilent Technologies, Inc.

## 概要

このアプリケーションノートでは、マレイン酸チモロールのキラル分離の開発と最適化について説明 します。このメソッドでは Agilent InfinityLab Poroshell 120 キラル-T カラムを使用し、マレイン酸 チモロールのキラル分離で用いられている現行の USP メソッドと比較します。

# はじめに

医薬品業界では、エナンチオマーの高速分離 がますます重要になっています。政府規制が強 化されるとともに、多数の光学活性医薬品化 合物が導入されているため、これらの化合物を 分析する際には、迅速かつ高感度で信頼性の 高いメソッドを考案することが重要になります。 現在使用されている薬物はその半分以上がキ ラル化合物であり、これらのほぼ90%が、2つ のエナンチオマーの等モル混合物で構成される ラセミ化合物として販売されています。キラル 薬物は化学構造は同じですが、その大部分 の異性体は生物学的活性が大きく異なってい ます。<sup>1,2</sup>

へキサンなどの順相溶媒による、セルロースま たはアミロースベースのキラル選択相(CSP) を用いた多数のキラル分離が実施されていま すが、メタノールなどのより一般的な溶媒を ベースにした分離では、他の相が要求される ことが多くなっています。これらの相は、逆相 メソッドを実施しているラボに簡単に組み込む ことができます。

InfinityLab Poroshell 120 キラル-T(テイコ プラニン)カラムは、逆相および順相 HPLC、 さらに SFC でも使用できます。また結合 CSP であるため、相のブリードや損失のリスクなし に、コーティング CSP 以外のさまざまな溶媒 で使用できます。

InfinityLab Poroshell 120 キラル-T およびそ の他のテイコプラニンベースの CSP は、極性 イオンモード(メタノールおよびアンモニウム 塩ベースの移動相、逆相(メタノールまたは アセトニトリルとバッファ))、または極性有機 モード(メタノール、エタノール、または別の 100 % 有機溶媒)で作用することがわかって います。逆相モードでは、有機成分および移 動相 pH の性質と濃度により、リテンションと 選択性が制御されています。<sup>4</sup> InfinityLab Poroshell 120 キラル-T (テイコ プラニン)のような糖ペプチドベースのキラ ルカラムは、逆相および順相 HPLC、さらに SFC のさまざまな溶媒で使用できます。糖ペ プチドは両性であり、イオン化した酸性および 塩基性グループの両方が含まれています。そ のため、糖ペプチドは移動相の pH に応じて、 正に帯電するか、負に帯電するか、または中 性になることができます。したがって、この種 の CSP を用いてイオン性化合物を分離する 際には、キラル認識にイオン性相互作用を関与 させることができます。この種の CSP における キラル認識では、今述べたことが主要な役割を 果たすと考えられています。キラル認識で CSP として抗生物質を使用することに関連して発生 する可能性があるその他の相互作用には、水 素結合、立体化、双極子-双極子、π-π相互 作用、および疎水性相互作用があります。こ れらの相互作用は、使用する個別の成分と移 動相モードの特性によって決定される、さまざ まな組み合わせにおいて発生する場合があり ます。各分離モードでは、キラル認識において 同時に異なる相互作用が発生します。これに より、多数のキラル分離およびこの種の CSP で適切に分離されるさまざまなタイプのキラル 化合物に関する情報が得られます。

糖ペプチドテイコプラニンは、表面多孔質シリ カ粒子と共有結合することにより、耐溶媒性 のある安定したクロマトグラフメディアを生成 します。この共有結合相は、一般的な HPLC 移動相およびメタノール、エタノール、IPA、 THF、リン酸塩、ギ酸塩、酢酸塩、ギ酸、TFA、 TEA、NH<sub>4</sub>OH などの添加物に対する耐性が あります。テイコプラニン結合相の構造を図 1 に示します。 このアプリケーションノートでは、InfinityLab Poroshell 120 キラル-T でのマレイン酸チモ ロールの分離を開発して最適化し、現行の USP メソッドと比較します。図2に、この化合 物の構造を示します。



**図 1.** Agilent InfinityLab Poroshell 120 キラル-T カラムでのテイコプラニン結合相の構造





# 実験方法

#### 試薬と実験方法

今回の実験では、Agilent 1260 Infinity II LC を使用しました。構成を表 1 に示します。基 本的なクロマトグラフィー条件を表 2 に示し ます。

マレイン酸チモロール、S-(-)-1-(t-ブチルアミ ノ)-3-[(4-モルホリノ-1,2,5-チアジアゾール-3-イル)オキシ]-2-プロパノールマレイン酸塩、欧 州薬局方 (EP) 参照標準、および R-チモロー ル欧州薬局方 (EP) 参照標準は、Sigma-Aldrich から購入しました。各エナンチオマー を 2 mg/mL でメタノールに溶解しました。 等しい量の各エナンチオマー溶液で構成され るサンプルを使用して、分離能を測定しまし た。LC/MS グレードのギ酸アンモニウムも Sigma-Aldrich から購入しました。HPLC グ レードのメタノールは Honeywell から購入し ました。2gのギ酸アンモニウムをメタノール に溶解し、0.2 % w/v で移動相を前処理しま した。低濃度の溶液は、この原液を希釈して 前処理しました。

# 結果と考察

4 本の InfinityLab Poroshell 120 キラルカラ ムと 6 つの移動相を用いたスクリーニングに より、チモロールキラル分離の初期分析を実 施しました。<sup>3,4</sup>このテストにより、0.2 % w/v ギ酸アンモニウム移動相含有のメタノールを 用いた 2 つのチモロールエナンチオマーとマ レイン酸塩の分離状態が明らかになりました。 この分離状態を図 3 に示します。

一般的に、極性イオン分離では、メタノールや アセトニトリルのような有機溶媒に少量の酸ま たは塩基を加えて使用します。または、今回の ケースのようにアンモニウム塩を使用できます。 ギ酸アンモニウムの濃度を低くすると、後半の 2 つのピークの分離能が高くなり、リテンション が増大することがわかりました。これらのピーク は、UV スペクトルを比較することにより、2 つの チモロールエナンチオマーとして同定しました。

#### 表1. 機器の構成

Agilent 1260 Infinity II LC			
Agilent 1260 Infinity II バイナリポンプ(G7112B)	最大 60 MPa、0 ~ 5 mL/min		
Agilent 1260 Infinity II マルチサンプラ(G7167A)	<ul> <li>・オートサンプラとヒーター:キャピラリー、ステンレス製、</li> <li>0.075 × 220 mm (p/n 5067-4784)</li> <li>・バイアル、スクリュートップ、茶色、ラベル付き、認定、2 mL、</li> <li>100 個 (p/n 5182-0716)</li> <li>・キャップ、スクリュー、青、PTFE/赤シリコンセプタム、</li> <li>100 個 (p/n 5182-0717)</li> </ul>		
Agilent Infinity II マルチカラムサーモスタット (G7116A)	<ul> <li>・超低分散ヒーター (G7116-60021)</li> <li>・ヒーターとカラム: Agilent InfinityLab クイックコネクトアセンブリ、 105 mm、0.075 mm (p/n5067-5961)</li> <li>・カラムと ELSD キャピラリー、ステンレス製、0.075 × 220 mm、 SV/SLV (p/n 5067-4784)</li> </ul>		
Agilent 1260 Infinity II DAD(G7115A)、 10 mm 光路長搭載	・G4212-6008、10 mm フローセル、1.0 μL V(ơ) ・40 Hz		
Agilent OpenLab CDS、バージョン C.01.07			

#### 表 2. LC メソッド条件

パラメータ	設定値	
カラム	Agilent InfinityLab Poroshell 120 $\pm \exists$ //-T, 4.6 $\times$ 100 mm, 2.7 $\mu m$ (p/n 685775-603)	
移動相	ギ酸アンモニウム含有メタノール (0~0.2%)	
流量	1 mL/min	
温度(カラム)	30 °C	
UV 波長	260 nm	
注入量	1μL	
サンプル濃度	2 mg/mL 水溶液	



図3.0.2% ギ酸アンモニウム含有メタノールを用いたマレイン酸チモロールの分離

この挙動は、極性イオン分離モードを用いた 際の代表的な分離です。イオン化グループが 存在しない分子のみをテイコプラニンまたは バンコマイシン相に加えると、メタノール、エ タノール、アセトニトリル、またはこれらの無 水溶媒の組み合わせで構成される移動相での 分離が高速化されます。極性有機分離の良い 例として、参考文献(キラルの概要)に示され ている 5-メチル-5-フェニルヒダントインまたは サリドマイドがあります。

ただし、化合物にイオン化グループが存在す る際には、少量の酸と塩基または揮発性塩を 加えることが必要になる場合があります。こ れは、大環状糖ペプチド自体の内部にイオン 化グループが存在するためです。後者の移動 相条件を極性イオンモードとして参照すること により、極性有機モードと区別しています。 極性イオンモードの選択性として、通常、化合物には最低2つの機能グループが存在しますが、そのうちの1つはイオン化されている必要があります。これらの機能グループには、アルコール、ハロゲン、いずれかの型の窒素(一次、二次、三次)、カルボニル、カルボキシル、酸化型の硫黄、またはリンを含めることができます。糖ペプチド相では、多数の酸と塩基、さらに揮発性塩を使用できます。酢酸アンモニウムは酸性分子の代表例であり、トリフルオロ酢酸アンモニウムおよびギ酸アンモニウム は塩基性分子で使用されます。チモロールのpKa は約9.2 であるため、MS 対応の塩であるギ酸アンモニウムの方が適しています。

最初のピークを、共通塩であるマレイン酸塩と して同定しました。図4に、ギ酸アンモニウム の3種類の濃度0.2、0.15、0.1%での分離 を示します。最初のチモロールのピークは、欧 州薬局方標準を注入することにより、「S」体 エナンチオマーとして同定しました。表3に、 6種類のギ酸アンモニウムの濃度、各濃度で の2つのチモロールエナンチオマーのリテン ション、およびチモロールエナンチオマーの分 離能を示します。これらのデータをプロットし

表 3. Agilent InfinityLab Poroshell 120 キラル-T でのさまざまなギ酸アンモニウム濃度におけるマレイ ン酸チモロールの分離

Rs	t1	t2
-	-	-
4.12	8.585	9.917
3.71	5.073	5.798
3.17	3.204	3.597
2.78	2.577	2.848
2.46	2.257	2.467
	<b>Rs</b> - 4.12 3.71 3.17 2.78 2.46	Rs         11           -         -           4.12         8.585           3.71         5.073           3.17         3.204           2.78         2.577           2.46         2.257



図4.0.2、0.15、0.1% ギ酸アンモニウム含有メタノールを用いたマレイン酸チモロールの分離

たものが図 5A と 5B です。図 6 に、この データセットのクロマトグラフの重ね表示を 示します。このデータは、ギ酸アンモニウムの 濃度が低くなると、チモロールエナンチオマー の分離能が高くなることを示しています。一方、 ピークのリテンションは増大しています。ギ酸 アンモニウムの濃度を低くすることにより、分 離能が高くなりました。ギ酸アンモニウムの濃 度 0.025 % w/v では、分離能は 4.12 です。 ギ酸アンモニウムの濃度 0 % では、カラムに より、チモロールのピークが識別できない場 所にピークが保持されています。

マレイン酸チモロール用の USP メソッドでは、 5 µm、4.6 × 250 mm 順相メソッドを利用し ており、USP 分類 L40 (セルローストリス-3,5-ジメチルフェニルカーバメートコーティング多孔 質シリカ) カラムとヘキサン/イソプロパノール/ ジメチルアミン、960:40:2 移動相が用いられ ています。このメソッドは、欧州薬局方にも記 載されています。この移動相にはジメチルア ミンが存在するため、MS には対応していま せん。<sup>5, 6, 7</sup>USP/EP メソッドに関するもう1つ の問題は、コーティングキラルカラムが使用さ れていることです。コーティングカラムは、固 定化キラルカラムと比較して寿命がかなり制 限されており、注入回数が 200 回程度の場合 もあります。アセトン、クロロホルム、DMF、 ジメチルスルホキシド、酢酸エチル、塩化メチ レン、THF など HPLC 溶出液で一般的に使 用されている多数の溶媒は、システム内の残 留物中に存在する場合でも、キラル固定相を 破壊することがあります。一般に、この順相溶 媒を使用しているシステムは、順相アプリケー ション専用です。<sup>8</sup>InfinityLab Poroshell 120 キラル-T は完全に結合されたカラムであり、 一般的に使用されているクロマトグラフィー溶 媒に対する耐性があります。

4.6 × 250 mm カラムで欧州薬局方メソッド を用いた分離について説明している参考文献 には、2つのチモロールエナンチオマーのリテン ションタイムが9~14分の間であることに加 えて、分離能4.8の要件に適合していること が示されています。同じカラムでSFCを用い ることにより、このメソッドを新しいメソッドと 比較しています。新しいメソッドは非常に高速 である一方、分離能が低下していることが報 告されています。<sup>9</sup>



図 5A. ギ酸アンモニウム濃度の関数としてのチモロールエナンチオマーのリテンションタイム



図 5B. ギ酸アンモニウム濃度の関数としてのチモロールエナンチオマーの分離能

### 結論

ここでは、極性イオン移動相(ギ酸アンモニウ ム含有メタノール)と InfinityLab Poroshell 120 キラル-T カラムを用いた、チモロールエ ナンチオマーの分離について説明しました。 今回のメソッドでは 220 nm UV 検出を用い ていますが、移動相は MS 検出と完全に互換 性があります。したがって、トラブルシューティ ングプロセスメソッドまたは法医学分析に簡 単に適合できます。このメソッドは、いずれの HPLC システムでも区別なく実施できます。



図 6. ギ酸アンモニウム濃度の関数としてのマレイン酸チモロールの分離

# 参考文献

- Bonner, W. A. Parity Violation and the Evolution of Biomolecular Homochirality. Chirality **2000**, 12, 114–126.
- Food and Drug Administration, FDA's Policy Statement for the Development of New Stereoisomeric Drugs. **1992**, 57 Fed. Reg. 22249.

### ホームページ

#### www.agilent.com/chem/jp

#### カストマコンタクトセンタ

## 0120-477-111 email\_japan@agilent.com

本製品は一般的な実験用途での使用を想定しており、 医薬品医療機器等法に基づく登録を行っておりません。 本文書に記載の情報、説明、製品仕様等は予告なしに 変更されることがあります。

アジレント・テクノロジー株式会社 © Agilent Technologies, Inc. 2020 Printed in Japan, June 23, 2020 5994-2143JAJP DE.0087615741

- Long, W. J.; Mack, A. E. Screening Chiral Compounds Using Superficially Porous Based Chiral Columns and a Method Development Solution. Pittcon 2018, CO-1479.
- 4. Put Agilent InfinityLab Poroshell 102 Chiral Innovations to Work on Your Challenging Separations. Agilent InfinityLab Poroshell 120 Chiral application compendium, publication number 5991-8450EN, **2017.**
- 5. United States Pharmacopeia USP 40 Timolol Maleate Monograph.
- 6. European Pharmacopoeia, 4th edition, Addendum 2002, Council of Europe, Strasbourg.

- 7 European Pharmacopoeia (Version 7.5): Timolol Maleate Monograph.
- Instruction Manual for Chiralcel OD-H and OJ-H Chiral Technologies 07/2013.
- Marley, A.; Connolly, D. Determination of (R)-Timolol in (S)-Timolol Maleate Active Pharmaceutical Ingredient: Validation of a New Supercritical Fluid Chromatography Method with an Established Normal Phase Liquid Chromatography Method. J. Chromatogr.A **2014**, 1325, 213– 220.

