

タンパク質凝集体の評価方法を最適化し 実行するための包括的なワークフロー

サイズ排除クロマトグラフィーと光散乱検出を組み合わせた メソッド開発

# アプリケーションノート

生物医薬品・バイオシミラー

## 著者

Andy Coffey and Matthew Rain Agilent Technologies, Inc

## 概要

このアプリケーションノートでは、包括的な凝集体分析のワークフローについて次の点を解説します。

- モノクローナル抗体の高性能なサイズ排除クロマトグラフィー (SEC) における移動相条件の 最適化
- 単量体、二量体、より高次の凝集体を含む凝集プロファイルの特性分析

今回の調査では、Agilent Buffer Advisor ソフトウェアを使用して、1260 Infinity II バイオイナート LC シ ステムのバイオイナートクォータナリポンプの機能をフル活用し、一連の高速 LC 分析中にリアルタ イムでさまざまな緩衝液の組成を自動的に調製し、複雑な SEC の最適化実験を自動化しました。 Agilent 1260 Infinity Bio-MDS マルチ検出器スイートはダイナミック光散乱検出機能を備え、高次のタ ンパク質凝集体を突き止め、分子量の絶対値を測定し、UV 検出システムによる定量測定を向上さ せます。



## はじめに

モノクローナル抗体 (mAbs) およびタンパク質 の中には溶液中で自然に凝集するものがあり ます<sup>12</sup>。多くのバイオ医薬品アプリケーション では、凝集の程度をさまざまな条件下で高い 精度で定性し定量する必要があります。サイ ズ排除クロマトグラフィー (SEC) は、タンパク 質凝集体の特性分析および定量のための優 れた手法ですが、正確に測定するには天然の タンパク質構造に対応する条件下での優れた クロマトグラフィーが必要です。特定のタンパ ク質の分離ピークの形状を改善し分離能を高 めるために、さまざまな移動相の条件を評価 することが多くの場合に必要となります。

緩衝液の条件の最適化の有用性は SEC 手法 で見落とされることがあります。従来より、緩 衝液の条件を最適化して固定相の材質との望 ましくない非特異的相互作用を克服していま したが、カラムに適した条件でないと、測定対 象とする凝集体そのものを攪乱するリスクを招 くことがあります。しかし、Agilent AdvanceBio SEC カラムの不活性表面コーティングは、さま ざまな緩衝液条件の二次的反応を抑制し、タ ンパク質の構造およびクロマトグラフィー分離 能が得られるように緩衝液の化学的性質を最 適化する優れた柔軟性を提供します。 凝集分 析のための SEC 緩衝液条件をルーチンで最適 化するうえでのもう一つの障害に、パラメータ スカウティング実験の多大な複雑さがありま す。最適化実験では移動相の可能性に関して 複雑な表の設計が必要となり、多数の異なる 溶液を手作業で作成し、塩、緩衝液、pHの変 化の組み合わせについて実験して評価しなけ ればなりません。しかし、最先端のツールによ り、SEC 条件を最適化するためのワークフロー を大幅に合理化して、凝集体の特性分析およ び定量分析を実行し、日常的に使用できる最 適化された手法を開発することができました。

このアプリケーションノートでは、凝集体の調 査における完全なワークフローソリューション の有用性を次の点について説明します。

- Agilent Buffer Advisor ソフトウェアと 高分解能バイオイナートクォータナリLC ポンプを用いた、4 種類のシンプルな 原液からの指定したリストのLC 緩衝液 の自動混合とpH および緩衝液濃度の リアルタイム調整
- ダイナミック光散乱検出による高次 タンパク質凝集体の測定、広い質量
   範囲にわたり感度を高めるための
   UV 検出での補完
- 光散乱検出による、凝集されたタンパク 質および単量体のタンパク質の分子量と 流体力学半径の絶対値の特性分析

今回の分析では、Agilent AdvanceBio SEC 15 cm カラムを使用して高速スクリーニングのため の高速分離を実現しました。AdvanceBio SEC カ ラムには、より長いものやより短いものなどさ まざまな長さのカラムが用意されているため、 追加の分離能が必要な場合には、30 cm のカ ラム、あるいは複数のカラムを組み合わせて 使用する場合にも容易にメソッドを適用できる ことが重要です。今回の調査で使用したモノ クローナル抗体サンプルには、市販のリツキ シマブおよびリツキシマブバイオシミラーが含 まれています。

## 実験方法

## 試薬、サンプル、材料

ー塩基性および二塩基性のリン酸水素ナトリ ウムと塩化ナトリウムは VWR から購入しまし た。化学物質および溶媒はすべて99.7 % 以 上の純度のものを使用しました。18 MΩ 未満 の水は、Milli-Q A10 純水生成装置 (Millipore、 米国) で生成しました。溶液は毎日調整し、 使用前に 0.22 μm メンブランフィルタでろ過し ました。

## 機器

凝集体分析には、次のモジュールで構成され る Agilent 1260 Infinity II バイオイナート LC シス テムを使用しました。

- Agilent 1260 Infinity II バイオイナート クォータナリポンプ (G5654A)
- Agilent 1260 Infinity II バイオイナート マルチサンプラ、サンプル冷却器および マルチウォッシュ搭載 (G5668A)
- Agilent 1260 Infinity II マルチカラム サーモスタット、バイオイナート 熱交換器搭載 (G7116A)
- Agilent 1260 Infinity II ダイオードアレイ 検出器 WR、バイオイナート標準 フローセル搭載 (G7115A)
- Agilent 1260 Infinity Bio-SEC マルチ検出器
  システム (G7805AA)
- Agilent AdvanceBio SEC 300Å、
  7.8 × 150 mm、2.7 μm カラム
  (p/n PL1180-3301) または
- Agilent AdvanceBio SEC、
  2.7 μm 300Å 7.8 × 300 mm カラム (p/n PL1180-5301)

### 機器条件

パラメータ	値
移動相	表1参照
流量	0.8 mL/min
温度	25 ° C
注入量	1 ~ 25 µL (サンプル濃度に依存)
検出	220 nm、280 nm、LS 90°、DLS
サンプル	リツキシマブ先発薬、
	リツキシマブバイオシミラー、BSA

## 結果と考察

各サンプル混合物についての最適な移動相の 組成を特定するために、4 つの pH レベルの 3 種類の移動相、つまり 12 の実験条件のマト リックスを作成し評価しました。緩衝液の組成 は次のとおりです。

- 150 mM リン酸ナトリウム
- ・ 10 mM リン酸ナトリウム + 140 mM NaCl (さまざまな pH 値での PBS のエミュレー ション)
- 100 mM リン酸ナトリウム + 150 mM NaCl

各移動相を、pH 6.2、6.6、7.0、7.4 でテストしま した。

各実験を行うために、Agilent Buffer Advisor で 各処理についての緩衝液の組成と pH を指定 しました。Agilent Buffer Advisor ソフトウェアに より、A ~ D 原液の適切な混合物を自動計算 し、LC パラメータスカウティング分析中に望ま しい移動相をリアルタイムで調製しました。表 1 に 12 の実験条件を示します。

これらの条件を使用してリツキシマブの市販 のサンプル、リツキシマブバイオシミラーのサ ンプル、市販の BSA タンパク質標準混合物 (機 器キャリブレーション用の 10 mg/mL 溶液) を 分析しました。

Agilent AdvanceBio SEC 300Å 150 × 7.8 mm カラ ムを使用することによって、1 サンプルあたり 10 分以内でスクリーニングできました。

#### 表1.実験の条件と対応する移動相の組成

		ユーザーが指定したパラメータ			ソフトウェアが計算したパラメータ			
		緩衝液	NaCl	合計濃度				
実験方法	pН	(mM)	(mM)	(mM)	% A	% B	% C	% D
1	6.2	150	0	150	25.0	0.0	57.0	18.0
2	6.6	150	0	150	25.0	0.0	42.3	32.7
3	7.0	150	0	150	25.0	0.0	26.3	48.7
4	7.4	150	0	150	25.0	0.0	13.8	61.2
5	7.4	10	140	150	67.0	28.0	0.9	4.1
6	7.0	10	140	150	67.0	28.0	1.8	3.2
7	6.6	10	140	150	67.0	28.0	2.9	2.1
8	6.2	10	140	150	67.0	28.0	3.9	1.1
9	6.2	100	150	250	20.0	30.0	36.3	13.7
10	6.6	100	150	250	20.0	30.0	26.1	23.9
11	7.0	100	150	250	20.0	30.0	15.6	34.4
12	7.4	100	150	250	20.0	30.0	7.8	42.2

A = 水

B = 500 mM NaCl

 $C=200 \text{ mM } \text{NaH}_2\text{PO}_4$ 

 $D = 200 \text{ mM} \text{ Na}_2\text{HPO}_4$ 

図1と図2に示すように、クロマトグラフィー データの最初のレビューから、実験中の移動 相の条件がリツキシマブ先発薬とリツキシマブ バイオシミラーのピークの形状に影響を与え ることが明らかになりました。2種類の分子は 似たプロファイルを示し、10 mM リン酸ナトリ ウムと140 mM NaCl の移動相組成では同様の 異常な挙動を示しています。この移動相の組 成は、ピーク高が低くなるとピークテーリング が著しく増大しました。このタイプの挙動は他 のタンパク質でも観察され、メソッド開発とメ ソッド堅牢性の両方を実現するために移動相 組成の影響を注意して評価する必要性を示し ています。



図 1.12 種類の異なる移動相条件で連続して、リツキシマブ先発品 (1 μL 注入) を分析 (表 1 の実験 1 ~ 12 を参照)。

凝集体の含有量の定量には、UV 検出器を使 用する必要があります。単量体のピークと凝 集体のピーク (検出された場所) を積分して、 ピーク面積の凝集体のパーセント率を求める ことができます。リツキシマブ先発薬の結果 を図 3A に、リツキシマブバイオシミラーの結 果を図 3B に示しています。150 mM リン酸ナ トリウムまたは 100 mM リン酸ナトリウムと 150 mM NaCl を用い pH 7.0 で、最も一貫性の ある結果が得られました (実験 3 と 11)。



図 2.12 種類の異なる移動相条件で連続して、リツキシマブバイオシミラー (1 µL 注入) を分析 (表 1 の実験 1 ~ 12 を参照)。

図 4 は、組成が異なる 3 種類の緩衝液で pH 7.0 で得られるクロマトグラムです。この分子 の分析に移動相として 10 mM リン酸ナトリウ ムと 140 mM NaCl を使用した場合には、凝集 体 (矢印部分) が検出されなかったことを明確 に示しています。

オプションで光散乱検出器を凝集体分析ワー クフローに追加すると、タンパク質凝集体につ いてより有益な情報が明らかになります。この BSA の場合、明確に定義された分子を1回注 入してシンプルな機器キャリブレーションを実 行した後、検出器間の遅延と機器定数を迅速 に求めることができます。Agilent Bio-SEC ソフト ウェアを用いると、同じ条件のカラムと流量で 得られた個別の任意クロマトグラムから個々 のピークについて分子量情報のレポートを作 成できます。



図 3A. リツキシマブ先発薬の凝集体含有量のピーク面積のパーセント率(表1の実験1~12を参照)。 凝集体が検出された実験のポイントが示されています。



図 3B. リツキシマブバイオシミラーの凝集体含有量のピーク面積のパーセント率(表1の実験1~12を参照)。 凝集体が検出された実験のポイントが示されています。



図 4. 最適化された pH 7.0 で異なる緩衝液塩濃度を用いたリツキシマブ先発薬 (A) と リツキシマブバイオシミラー (B) 分析の UV 220 nm 信号のベースラインの拡大図 (表 1 の実験 3、6、11 を参照)。

図5は、実験11の高塩濃度条件下でのリツ キシマブ先発薬とリツキシマブバイオシミラー の光散乱分析の結果を示しています。単量体 のピークのみの分析に焦点を当てると、光散 乱分析により以前のアプリケーションノート<sup>1</sup> に記載されている精密質量値に近い分子量が 報告されました。バイオシミラー分子は、先発 薬分子でははっきりと見られないC末端リジ ン変異体が存在しているため、質量がわずか に大きくなることが予想されます。

より大きな凝集体と肉眼では見えない粒子の 形成は、バイオ医薬品アプリケーションで特に 懸念していることです。高度に凝集されたサン プルに対する光散乱検出の感度は、UV 検出か ら得られた濃度データを補完します。

UV 検出による単量体と二量体の分析結果は きわめて類似しています。しかし、図6が示す ように、特定の移動相の条件下では光散乱検 出器はより高次の凝集体により大きく反応す ることが立証され、リツキシマプ先発薬とバイ オシミラーのより広い範囲の凝集体ではいくつ かの相違点があることが明らかになりました。



図 5.100 mM リン酸ナトリウムと 150 mM NaCl、pH 7.0 でのリツキシマブ先発薬 (A) と リツキシマブバイオシミラー (B) の LS 分析の比較 (表 1 の実験 11)。



図 6. 150 mM リン酸ナトリウム、pH 7.0 でのリツキシマブ先発薬 (A) とリツキシマブバイオシミラー (B) の LS 分析の比較 (表 1 の実験 3)。

DLS 機能の追加は得られる情報の内容をさら に詳細にして、流体力学半径の測定を可能に します (図 7)。

## 結論

完全にバイオイナートな流路の Agilent 1260 Infinity II Bio-inert LC システムと Agilent Buffer Advisor ソフトウェアを組み合わせると、タンパ ク質凝集体の定量のためのサイズ排除クロマ トグラフィーに適したメソッドの最適化を簡単 に実現できます。より短い Agilent AdvanceBio SEC 300Å 15 cm カラムは分離を高速化し、ス ループットを大幅に高め広い範囲の分析条 件のスクリーニングに必要な時間を短縮しま す。分離能と真度を高めるには、より長い30 cm カラムを使用します。AdvanceBio SEC カラ ムには、非特異的な結合がほとんど生じない などの利点があります。Bio-MDS マルチ検出 器スイートと Bio-SEC ソフトウェアを使用すれ ば、他のどの手段でも検出が困難な低濃度の 高分子量凝集体を突き止めることができます。 最終的に、Bio-MDS を使用して、タンパク質分 子量を測定したり、DLS 検出と組み合わせて 流体力学半径情報を求めることができます。 いくつかの技術の組み合わせにより包括的な ワークフローソリューションを構成して、SEC 条件を短時間で最適化し、適切な緩衝液条件 で凝集体を分子量範囲全体で正確に定量し、 モノクローナル抗体の凝集動態の特性評価を 行うことができます。



図 7.LS 検出器による流体力学半径の結果を示すリツキシマブバイオシミラー DLS 分析。

## 参考文献

- Guidance for Industry Immunogenicity Assessment for Therapeutic Protein Products, U.S. Department of Health and Human Services Food and Drug Administration Center for Drug Evaluation and Research (CDER) Center for Biologics Evaluation and Research (CBER), 2014.
- Mahler, H-C.; et al. Protein Aggregation: Pathways, Induction Factors and Analysis, J. Pharm. Sci. 2008, 98(9).
- Schneider, S. 2D-LC/MS Characterization of Charge Variants Using Ion Exchange and Reversed-Phase Chromatography, *Agilent Technologies Application Note* publication number 5991-6673EN 2016.

#### 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。アジレントは、本文書に 誤りが発見された場合、また、本文書の使用により 付随的または間接的に生じる損害について一切免責と させていただきます。

アジレント・テクノロジー株式会社 © Agilent Technologies, Inc. 2017 Printed in Japan, January 5, 2017 5991-7476JAJP

