

モノクローナル抗体の複数の 重要品質特性評価および 比較可能性評価

著者

Suresh Babu C.V., Brian Liau Agilent Technologies, Inc.

概要

多特性解析メソッド (Multi-attribute method:MAM) に基づくLC/MS ペプチドマッピングは、バイオ 医薬品分析の分野で大きな注目を集めています。本アプリケーションノートでは、さまざまなモノク ローナル抗体 (mAb)の重要品質特性 (CQA)を特定・定量するためのLC/MS ペプチドマッピング アプローチについて説明します。このペプチドマッピングメソッドを従来の方法と比較したところ、リジン 切断、酸化、脱アミド化、N 末端環化、およびグリコシル化に関して、非常に類似した推定値が得ら れることがわかりました。

はじめに

モノクローナル抗体 (mAb) は生体系に由来 するもので、急速な成長を遂げている抗体分子 です。この複雑な分子は生物由来であるため、 生産中にさまざまな翻訳後修飾(PTM)を受 け、不均一性がかなり高くなります。 mAb の開 発および製造中に、これらの意図しない変化を CQA として特定し、監視することが重要です。 この目的のために、各 CQA の化学的な相違を 利用した分析試験方法がいくつか伝統的に用 いられていますが、各技法にはさまざまな長所 と短所があります。最近では、LC/MS ベース のマルチ特性メソッド (MAM) ¹ アプローチが 広く用いられるようになってきています。 mAb CQA を包括的に特定する単一のメソッドとし て利用することができ、従来のメソッドと比べ て分析時間とコストを大幅に節約できるから です。ただし、規制の厳しい品質管理環境に MAM アプローチを適用する前に、まず、従来 の分析方法がペプチドマッピングの結果と正確 に一致することを確認する必要があります。ペ プチドマッピングのワークフローにより、従来の アッセイと同等の結果が得られることを確認す ることが重要です。

本研究では、ペプチドマッピングメソッドの性 能を、イオン交換クロマトグラフィー、親水性 相互作用クロマトグラフィー(HILIC)、サブユ ニットレベルの逆相クロマトグラフィーなどの 従来のアッセイと比較します。mAbのペプチ ドマッピングは、Agilent AssayMAP Bravo Liquid Handling Platform、Agilent 1290 Infinity II LC システム、Agilent 6545XT AdvanceBio LC/Q-TOF、および Agilent MassHunter BioConfirm ソフトウェアによ るデータ解析から構成される統合型ワークフ ローを使用しました。各 CQA の相対 % 定量 結果を、ペプチドマッピングメソッドと従来の アッセイとの間で比較しました。

装置、カラム、ソフトウェア

サンプル前処理自動化システム Agilent AssavMAP Bravo Liquid Handling Platform (G5542A) LC システム Agilent 1260 Infinity II バイオイナート LC Agilent 1290 Infinity II LC システム ・Agilent 1260 Infinity II バイオイナートポンプ (G5654A) ・Agilent 1290 Infinity II ハイスピードポンプ (G7120A) ・ Agilent 1260 Infinity II バイオイナートマルチサンプラ • Agilent 1290 Infinity II マルチカラムサーモスタット (G5668A) サンプル冷却器付 (G7116B) • Agilent 1260 Infinity II マルチカラムサーモスタット ・Agilent 1290 Infinity II マルチサンプラ (G7167B) (G7116A) バイオイナート熱交換器付 ・Agilent 1260 Infinity バイオイナート ・分析スケールフラクションコレクタ(G5664A) 検出 ・Agilent 1260 Infinity II ダイオードアレイ検出器 WR (G7115A) バイオイナートフローセル付 · Agilent 1260 Infinity 蛍光検出器(G1321B) Agilent 6545XT AdvanceBio LC/Q-TOF カラム ・Agilent Bio MAb、非ポーラス、4.6 imes 250 mm、5 μ m HPLC、PEEK Agilent ZORBAX RRHD 300 SB-C18, 2.1 × 100 mm, 1.8 um ・Agilent AdvanceBio グリカンマッピング、2.1 × 150 mm、1.8 μm ・Agilent AdvanceBio ペプチドマッピング、2.1 imes 150 mm、2.7 μ m、120 Å ソフトウェア ・Agilent OpenLab CDS バージョン 2.3 Agilent Buffer Advisor A.01.01 [009] ・LC システム用 Agilent OpenLAB CDS ChemStation Edition、Rev. C.01.07 Agilent MassHunter Acquisition (B.10.00) • Agilent MassHunter BioConfirm 10.0

Agilent VWorks Automation Control 13.1.1.1383

試料調製

mAb1、mAb2、および mAb3 は地元の販売 代理店(シンガポール)から購入し、製造元 の指示に従って保管しました。トリズマ塩基、 トリス-HCl、グアニジン塩酸塩、トリス(2 カ ルボキシエチル)ホスフィン(TCEP)、ヨード アセトアミド(IAA)、ギ酸、トリフルオロ酢酸、 リン酸ーナトリウム、リン酸ニナトリウム、塩 化ナトリウム、および LC/MS グレードの溶媒 は、Sigma-Aldrich から購入しました。IdeS プロテアーゼは Genovis から購入しました。 高品質配列グレードのトリプシンは、Agilent Technologies で用意しました。

実験メソッド

電荷変異体: サンプルは希釈せず直接注入しま した(10 mg/mL)。表1 および2 に、mAb1 および mAb2 の弱カチオン交換(WCX) クロ マトグラフィーで使用したクロマトグラフィーパ ラメーターを示します。収集した WCX フラク ションに対して、トリプシン消化およびペプチド マッピング分析を行いました。

酸化:5 µL の mAb サンプル (2 mg/mLト リス-HCl バッファ)を5 µL の ldeS プロテ アーゼに添加し、30 ℃で 30 分間インキュ ベートした後、分析の前に室温まで放冷しま した。 **遊離グリカン:** 標識 N-グリカンサンプルは、 GlykoPrep 高速 N-グリカンキット(GPPNG-PC)を用いて、AssayMAP Bravo Liquid Handling システムを使用して調製しました。² 標識 N-グリカンは HILIC で分離しました。

トリプシン消化

フラクションを収集した WCX サンプルと mAb サンプルを還元/アルキル化し、トリプシ ンで消化した後、Agilent AssayMAP Bravo Liquid Handling Platform を用いて脱塩しま した。³ mAb を含むサンプルプレートを変性 バッファ(8 M グアニジン-HCl、5 mM TCEP と150 mM トリス pH8)で再構成し、60°C で 60 分間インキュベートしました。ジスルフィド 結合を変性および還元した後、133 mM ヨード アセトアミドを添加して遊離システインのアル キル化を行いました(室温で 40 分)。その後、 1.75 % TFA でサンプルを酸性化し希釈しまし た。さらに、サンプルを RP-W カートリッジを 使用して Protein Cleanup アプリケーション にかけ、37 °C で一晩トリプシン消化(20:1、 タンパク質:プロテアーゼ w:w)を行いまし た。その後、AssayMAP 試薬トランスファー ユーティリティを用いてサンプルを 0.1 % ギ酸 で酸性化してトリプシン活性を停止しました。 AssayMAP Bravo Peptide Cleanup プロト コル(脱塩)は、C18 逆相カートリッジを使 用しました。

表 1. LC の分析条件

	WCX (C 末端リジン、N 末端環化変異体、 脱アミド化)	RP (酸化)	HILIC (遊離グリカン)	RP (ペプチドマッピング)
パラメータ	Agilent 1260 Infinity II バイオイナート LC	Agilent 1290 Infinity II LC	Agilent 1290 Infinity II LC	Agilent 1290 Infinity II LC
カラム	Agilent Bio MAb NP5、 4.6 \times 250 mm、5 μ m PEEK	Agilent ZORBAX RRHD 300 SB-C18、 2.1 × 100 mm、1.8 μm	Agilent AdvanceBio グリカンマッピング、 2.1 × 150 mm、1.8 μm	Agilent AdvanceBio ペプチドマッピング、 2.1 × 150 mm、2.7 µm、120 Å
注入量	$1\sim 10\mu L$	4 µL	2 µL	2 ~ 8 µL
サンプルの サーモスタット	5 ℃	5 ℃	5 °C	5 ℃
移動相	A) 水 B) NaCl (1,000 mM) C) NaH₂PO₄ (55 mM) D) Na₂HPO₄ (50 mM)	A) 0.1 % FA、0.1 % TFA 水溶液 B) IPA 80 %、ACN 10 %、水 10 % の 混合溶媒中 0.1 % FA、0.1 % TFA	A) 50 mM ギ酸アンモニウム、 pH 4.5 B) ACN	A) 0.1 % ギ酸水溶液 B) 0.1 % ギ酸 ACN 溶液
グラジエント	mAb1 まよじまととととととのでした。 時間(分) %A %	0 分 → 20 % B 2 分 → 20 % B 42 分 → 35 % B 42.5 分 → 80 % B 45 分 → 80 % B 45.1 分 → 20 % B	0 分 → 25 % B 2 分 → 30 % B 15.5 分 → 37 % B 17 分 → 80 % B 20 分 → 80 % B 21 分 → 20 % B 30 分 → 25 % B	0 分 \rightarrow 3 % B 30 分 \rightarrow 40 % B 33 分 \rightarrow 90 % B 35 分 \rightarrow 90 % B 37 分 \rightarrow 3 % B 40 分 \rightarrow 3 % B
カラム温度	25 °C	74 °C	40 °C	60 °C
流量	 0.75 mL/分	0.4 mL/分	15.5 ~ 20 分0.25 mL/分、それ以外は 0.4 mL/分	0.4 mL/分
UV	220/280 nm	-		-
FLD	-	-	Ex/Em = 285 nm/345 nm	-
フラクション コレクションモード	時間ベースのフラクショントリガ	-	-	-

表 2. LC/MS の条件

	WCX (C 末端リジン、						
パラメータ	N 末端環化変異体、脱アミド化)	RP(酸化)	HILIC (遊離グリカン)	RP (ペプチドマッピング)			
Agilent 6545XT AdvanceBio LC/Q-TOF							
イオンモード	-	ポジティブイオンモード、デュアル AJS ESI	ポジティブイオンモード、デュアル AJS ESI	ポジティブイオンモード、デュアル AJS ESI			
ドライガス温度	-	350 °C	150 °C	325 ℃			
ドライガス流量	-	12 L/分	9 L/分	13 L/分			
シースガス温度	-	350 °C	300 °C	275 ℃			
シースガス流量	-	11 L/分	10 L/分	12 L/分			
ネブライザ	-	35 psi	35 psi	35 psi			
キャピラリー電圧	-	4,000 V	3,000 V	4,000 V			
ノズル電圧	-	2,000 V	500 V	0 V			
フラグメンタ電圧	-	180 V	180 V	175 V			
スキマー電圧	-	65 V	65 V	65 V			
リファレンス質量	-	922.009798	922.009798	121.050873、922.009798			
取り込みモード	-	拡張(10,000 m/z)質量範囲	低質量、高分解能(4 Ghz)	拡張ダイナミックレンジ (2 GHz)			
MS 質量範囲	-	800~5,000 <i>m/z</i>	300~1,700 <i>m/z</i>	110~1,700 <i>m/z</i>			
取り込みレート	-	1 スペクトル/秒	2 スペクトル/秒	8 スペクトル/秒			
自動 MS/MS 範囲	-	-	-	50~1,700 <i>m/z</i>			
MS/MS 取り込みレート	-	-	-	3 スペクトル/秒			
選択幅	-	-	-	狭い(~ 1.3 <i>m/z</i>)			
プリカーサ/サイクル	-	-	-	上位 10			
コリジョンエネルギー	-	-	-	 荷電状態 傾き オフセット 2 3.1 1 3 および >3 3.6 -4.8 			
MS/MS の スレッシュホールド	-	-	-	1,000 カウントと 0.001 %			
ダイナミック排除オン	-	-	-	1 回繰り返し、0.1 または 0.2 分間排除			
プリカーサアバンダンス ベースのスキャンスピード	-	-	-	使用する			
目標値	-	-	-	25,000 カウント/スペクトル			
MS/MS 累積時間制限 の使用	-	-	-	使用する			
純度	-	-	-	100 % ストリンジェンシー、30 % カットオフ			
同位体モデル	-	-	-	ペプチド			
プリカーサの順位	-	-	-	荷電状態の次に存在量: +2、+3、>+3			

結果と考察

C 末端リジンの切断

■ 重鎖の C 末端リジンの切断は、カルボキシ ペプチダーゼ活性により一般的に観察される PTM です。リジン残基が失われると、mAb の正味の正電荷が減少し、電荷が不均一にな ります。C 末端リジン塩基性変異体(Lys あり、 Lys なし) は、カチオン交換クロマトグラフィー (CEX) でうまく分離できます。図 1A は、 Agilent Bio MAb PEEK カラムで高分解能分 離した mAb1 の電荷変異体プロファイルを示し ています。3つの異なるピークがあることがわか ります。13.11 分のピークをメインピーク(M) とします。早い溶出ピークと遅い溶出ピーク を、それぞれ酸性変異体(A1、A2)、塩基性 変異体 (B1、B2、B3、B4、B5、B6) とします。 メイン、酸性、および塩基性の電荷変異体の ピーク面積パーセントを図 1A 中の表に示し ます。LC/MS ペプチドマッピング分析によっ て C 末端リジンの切断を特性解析する前に、 電荷変異体ピークを個別のフラクションとして収 集し、複数回の注入にわたってプールしました。 ペプチドマッピング分析に基づいて、CEX ピー クを C 末端リジン切断変異体でアノテートし ます。C 末端リジンの不均一性のさらなる確 認を、カルボキシペプチダーゼ B (CPB) 処理 によって行いました。C 末端リジン切断の相対 含有量は、電荷変異体ピークの面積パーセント から計算し、58.28%であることがわかりました (表3)。

C 末端リジン変異体プロファイルのペプチド マッピングベースの定量は、mAb1をLC/MS ペプチドマッピングメソッドで分析すること で行いました。図 1B に mAb1 の LC/MS ペプチドマッピング分析の結果を示します。 MS/MS 分析により、修飾および未修飾の重 鎖 C 末端リジン切断ペプチドを同定しました。 相対含有量は 55.30 % であることがわかりま した。 従来の CEX およびペプチドマッピングメソッド の C 末端リジン切断変異体により定量した結 果の比較を表3に示します。どちらの方法でも、 C 末端のリジン切断の相対量の推定値は類似 したものでした。







図 1B. 2.1 × 150 mm、2.7 µm カラムを使用した mAb1 の Agilent LC/MS ペプチドマッピング

表 3. 相対 % 定量比較(従来のマッピング手法と ペプチドマッピングメソッド)

手法	相対含有量 %					
C 末端リジン変異体						
CEX	58.28					
ペプチドマッピング	55.30					
N 末端環化変異体(HC)						
CEX	100					
ペプチドマッピング	99.87					
Asp55 脱アミド化						
CEX	3.86					
ペプチドマッピング	4.79					
酸化(M256 + M432)						
RP-LC/MS (サブユニット)	2.38					
ペプチドマッピング	2.33					
グリカン						
	HILIC	ペプチドマッピング				
Man5	10.9	12.43				
G0	6.48	8.35				
G0F	49.13	50.41				
G1F	28.80	24.57				
G2F	4.69	4.24				

N 末端環化

N 末端の環化は、N 末端の Glu または Gln がピログルタミン酸 (pyro-Glu) に転位する ことによって引き起こされます。これにより 正味の正電荷が減少し、mAb はより酸性に なります。C 末端リジンの切断と同様に、こ の PTM に起因する電荷の不均一性は、CEX により確認できます。 図 2A に、Agilent Bio MAb PEEK カラムで mAb2 の電荷変異体を 高分解能分離した結果を示します。13.1 分の ピークをメインピーク(M)とします。早い溶 出ピークと遅い溶出ピークを、それぞれ酸性 変異体(A1、A2)、塩基性変異体(B1、B2、 B3) とします。 図 2A は、メイン、酸性、およ び塩基性電荷変異体の面積パーセントをまと めたものです。LC/MS ペプチドマッピング分 析を行って N 末端環化変異体を特性解析する 前に、電荷変異体ピークを個別のフラクション として収集し、複数の注入にわたってプールし

ました。ペプチドマッピング分析に基づいて、 CEX ピークを N 末端環化変異体でアノテート します。重鎖 N 末端環化の相対含有量は、電 荷変異体ピークの面積パーセントから計算し、 100%であることがわかりました(表 3)。

重鎖 N 末端環化のペプチドマッピングベース の定量は、mAb2 を LC/MS ペプチドマッピン グメソッドで分析することで行いました。図 2B に mAb2 の LC/MS ペプチドマッピング分析 の結果を示します。MS/MS 分析により、重鎖 ペプチドの N 末端環化を確認し、N 末端環化 を示す重鎖分子の割合が 99.98 % であること がわかりました。

従来の CEX とペプチドマッピングメソッドの N 末端環化変異体による定量の比較を表 3 に示します。どちらの方法でも、N 末端環化 の相対量の推定値は類似したものでした。



図 **2A.** Agilent 1260 Infinity バイオイナートクォータナリ LC で測定した mAb2 の電荷変異体プロファイル。 Agilent Bio Mab、4.6 × 250 mm、5 µm、PEEK カラムを使用



図 2B. 2.1 × 150 mm、2.7 µm カラムを使用した mAb2 の Agilent LC/MS ペプチドマッピング

脱アミド化

脱アミド化は mAb で最も頻繁に起こる PTM であり、タンパク質分解経路に関連しています。 アスパラギン (Asn、N)残基の脱アミド化で は、Asn が酸性異性体アスパラギン酸 (asp、 D)とイソアスパラギン酸 (isoAsp)に変換さ れ、正味の正電荷が減少します。CDR 領域の Asn 脱アミド化により、酸性の電荷変異体が 生じます。図 3A に、Agilent Bio MAb PEEK カラムで mAb3 を高分解能分離した際の電 荷変異体プロファイルを示します。11.4 分の ピークをメインピーク (M)とします。早い溶 出ピークと遅い溶出ピークを、それぞれ酸性 変異体 (A1、A2)、塩基性変異体 (B1、B2、 B3、B4)とします。図 3A は、メイン、酸性、 および塩基性電荷変異体の面積パーセントを まとめたものです。LC/MS ペプチドマッピング 分析を行って脱アミド化を特性解析する前に、 電荷変異体ピークを個別のフラクションとして 収集し、複数の注入にわたってプールしました。 ペプチドマッピング分析により、フラクション A2 の Asn55 脱アミド化を同定しました。N55 脱 アミド化の相対含有量は、電荷変異体ピーク の面積パーセントから計算し、3.86 % である ことがわかりました(表 3)。 脱アミド化のペプチドマッピングベースの定 量は、mAb3 を LC/MS ペプチドマッピング メソッドで分析しました。図 3B に mAb3 の LC/MS ペプチドマッピング分析の結果を示 します。MS/MS 分析により、修飾および未 修飾の IYPTN⁵⁵GYTR ペプチドが同定され、 Asn55 脱アミド化の相対含有量は4.79 % で あることがわかりました。

従来の CEX およびペプチドマッピングメソッド による脱アミド化の定量の比較を表 3 に示し ます。どちらの方法でも、Asn55 脱アミド化レ ベルの相対量の推定値は類似したものでした。



図 3A. Agilent 1260 Infinity バイオイナートクォータナリ LC で測定した mAb3 の電荷変異体プロファイル。 Agilent Bio Mab、4.6 × 250 mm、5 µm、PEEK カラムを使用



図 3B. 2.1 × 150 mm、2.7 µm カラムを使用した mAb3 の Agilent LC/MS ペプチドマッピング

酸化

メチオニン (Met) の酸化は mAb では一般 的であり、機能変化の可能性があるタンパク 質の構造変化をもたらします。Fc 領域には、 2 つのメチオニン (Met) 残基 (M256 およ び M432) が含まれています。これらは酸化 を受けやすい部位です。酸化は +16 Da の質 量増加をもたらし、ペプチドとインタクト mAb の両方で測定できます。

図 4A は、Agilent ZORBAX RRHD 300 SB-C18 カラムを使用して分離された Ides 消化 mAb2 Fc サブユニットの LC/MS プロファイ ルを示しています。デコンボリュートしたした 質量スペクトルの主要なピークは、2 つの酸 化種に対応します。これらは主に M256 と M432 に起因するものです。Fc 酸化の定量レ ベルは、デコンボリュートした質量スペクトル のピーク面積から計算し、2.38 % であること がわかりました。

酸化のペプチドマッピングベースの定量は、 mAb2 を LC/MS ペプチドマッピングメソッドで 分析することで行いました。 図 4B に、トリプシン mAb2 ペプチドの LC/MS プロファイルを示 します。MS/MS 分析により、酸化および未 修飾のペプチド変異体 (DTLM²⁵⁶ISR および WQQGNVFSCSVM⁴³²HEALHNHYTQK) が 同定され、相対的な酸化量 (M256 + M432) は 2.33 % であることがわかりました。

表3に、サブユニット分析またはペプチドマッ ピングメソッドのいずれかを使用して酸化の定 量を行い比較した結果を示します。どちらの 方法でも同様の推定値が得られました。



図 **4A.** Agilent ZORBAX RRHD 300 SB-C18、2.1 × 100 mm、1.8 μm、PEEK カラムを使用した Ide 消化 mAb2 のLC/MS 分析。Fc フラグメントのデコンボリュートした質量スペクトルは、32 Da のシフトを示します (2 つの酸素添加に相当)。





グリコシル化

アスパラギン結合 N 結合グリコシル化は、最も 重要な PTM の 1 つで、mAb の不均一性に 寄与します。グリカンの特性解析は、バイオ医 薬品の製造において極めて重要です。図 5A に、 Agilent AdvanceBio グリカンマッピングカラム で分離した mAb1 の InstantPC 標識 N-グリ カンの蛍光クロマトグラムを示します。典型的 な主要グリカン GOF、G1F、および G2F が示 されています。グリカン種の分離は、相対含有 量を計算するのに十分なものです。挿入表に、 各グリカンの相対量を示しました。グリカンの アノテーションを確認するために、各ピークか らの MS シグナルを、多価 H⁺ および Na⁺ 付 加イオンの同位体モデルと照合しました。

グリカン種のペプチドマッピングベースの 定量のために、mAb1 のトリプシンペプチ ドを LC/MS で分析しました。図 5B は、 AdvanceBio ペプチドマッピングカラムを使 用して分離されたさまざまなグリコシル化 TKPREEQYNSTYR ペプチドの抽出イオンク ロマトグラム (EIC)の重ね合わせ図を示して います。MS 分析により、TKPREEQYNSTYR ペプチドの主要なグリコフォーム (G0、G1F、 G0F、および G2F)と一致する質量が同定 されました。最も量の多いグリコフォームは、 G0F および G1F に対応します。

遊離グリカン (HILIC) とペプチドマッピング メソッドによるグリカン定量の比較を図 5C と 表 3 に示します。どちらの方法でも同様の推 定値が得られました。相関計数は 0.98 であ り、主要なグリカンに関しては、HLIC と RP ペプチドマッピングメソッドの間には強い正の 相関があることがわかります。











図 5C. HILIC 法とペプチドマッピングメソッドによるグリカン定量の比較をした相関グラフ

結論

本研究では、mAb の複数の CQA をモニタ リングするための Agilent ペプチドマッピング ワークフローについて説明しまた。ペプチド マッピングと従来のアッセイの間で複数の特 性の定量値の相対%値を比較したところ、非 常に良好な相関関係が示されました。結論と して、ペプチドマッピングアプローチは、従来 のアッセイの代替となり、mAb の品質特性の モニタリングに用いることができます。これに より、分析時間の短縮とコスト削減が可能 です。

参考文献

- Rogers, R. S. et al. A View on the Importance of "Multi-Attribute Method" for Measuring Purity of Biopharmaceuticals and Improving Overall Control Strategy. *The AAPS Journal* 2018, 20, 7.
- 2. A Comparative Study of the Intact Mass, Subunit Mass, and Released Glycans of Two Rituximab Biosimilars Using High-Resolution LC/MS, *Agilent Technologies application note*, publication number 5994-1653EN, **2020**.
- Rapid Antibody Digestion Enabled by Automated Reversed-Phase Desalting on the Agilent AssayMAP Bravo Platform, *Agilent Technologies application note*, publication number 5991-6478EN, **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・テクノロジー株式会社 © Agilent Technologies, Inc. 2020 Printed in Japan, August 19, 2020 5994-2245JAJP DE.0531828704

