バイオ医薬/医薬



シングル四重極 LC/MS を使用した、 インタクトおよびサブユニットレベルでの 抗体のグリコシル化のモニタリング

著者

Stephen Sciuto Agilent Technologies, Inc.

概要

このアプリケーションノートでは、Agilent InfinityLab Pro iQ Plus シングル四重極質量分析計(MS)を 用いた、トラスツズマブのインタクトおよびサブユニットレベルにおける抗体のグリコシル化の相対アバン ダンスのモニタリングについて説明します。インタクトレベルでは、アジレントの高分解能精密質量シス テムで測定済みのものと一致する相対アバンダンスにより、MS スペクトルデコンボリューションソフト ウェアから 5 種類のグリコシル化成分を同定しました。サブユニットレベルでは、トラスツズマブの軽 鎖はアミノ酸配列に基づき予想される単一成分を示し、一方でトラスツズマブの重鎖により G0、G0F、 G1F、G2F グリカンと一致する 4 つのピークが同定されました。グリコシル化モニタリングのための 信頼性の高い相対アバンダンスと、優れた MS スペクトルピーク形状の両方が、Agilent Pro iQ Plus 質量分析システムの性能を際立たせています。

はじめに

グリコシル化は、バイオ医薬品分野において モニタリングすべき重要なパラメータです。グリ カンは、医薬品の構造安定性、安全性、効能 に影響を及ぼすからです。¹許容範囲に収まる タンパク質グリコシル化の相対アバンダンスを 確保することが一般的な品質指標であり、繰り 返しモニタリングされます。相対アバンダンスが 品質管理(QC)チームがあらかじめ規定した 標準に合格しない場合、医薬品の製造プロセス に対してさらなる調査が必要になります。質量 分析は、モノクローナル抗体(mAb)の特性 解析のための感度と堅牢性に優れたメソッド を提供するため、グリカンの特性解析に最適 です。

この研究では、Agilent InfinityLab Pro iQ Plus 液体クロマトグラフィー /質量分析法(LC/ MS)と、優れた質量範囲および質量スペク トルピーク形状を使用し、インタクト/サブユ ニットレベルにおけるグリコシル化を検出しま した(HC、LC)。デコンボリューション後の グリコシル化の相対アバンダンスは、この機器 プラットフォームでタンパク質グリコシル化に 関する重要な情報²を評価できることを示して います。

実験方法

化学物質と溶液の調製

還元反応のために、Sigma-Aldrich(米国ミ ズーリ州セントルイス)の以下の溶液を使用し ました。

- A. 8 M 塩酸グアニジン、pH 8.5 (部品番号 G7294、100 mL)
- B. 0.2 M DTT (DL-ジチオスレイトール、部品番号 D9779、1 g)
- C. 1 M トリスバッファ、pH 8.0 (部品番号 648314, 100 mL)

変性:6.4 M グアニジン-HCl、200 mM トリス -HCl バッファ、pH 8.1

100 mL 計量フラスコで 80 mL の 8.0 M グア ニジン HCI 溶液を 20 mL の 1 M トリス-HCI バッファに添加します。反転により完全に混 合してから、ピペットと pH 試験紙を使用して pH を測定し、記録します。pH が 7.2 ~ 8.5 の 範囲に収まる場合は、還元ステップに進んで かまいません。

還元:50 mM トリス-HCl 溶液中の 200 mM DTT

分析用天秤を使用して 31 mg の DTT を測定 し、1 mL の 50 mM トリス-HCI 溶液を添加し て溶解します。還元ステップの前に溶液をボル テックスします。

標準とサンプル前処理

22 mg/mL の濃度のモノクローナル抗体 (mAb) トラスツズマブは Genentech (カリ フォルニア州サウスサンフランシスコ)から調 達しました。210 μ L の 0.1 % ギ酸水溶液に 10 μ L の原液を添加することにより、この溶液 を 1 μ g/ μ L の濃度に希釈しました。55 μ L の 1 μ g/ μ L 溶液を添加して、さらに 165 μ L の 0.1 % ギ酸水溶液を添加することにより、LC/ MS 分析に使用する最終溶液を 250 ng/ μ L の 濃度に希釈しました。

トラスツズマブの還元に使用したサンプル前処 理の手順:

- 2.3 µLのトラスツズマブ原液(22 mg/mL) を 27.4 µLの 6.4 M 塩酸グアニジン、 200 mMトリス-HCI に添加します。
- 50 mM トリス-HCI 中の 6.3 µL の
 200 mM DTT をトラスツズマブ溶液に添加します。静かに混ぜてスピンダウンしてから、37 ℃ で 30 分間(または60 ℃ で30 分間)インキュベートします。
- 3. 還元反応を停止するために、溶液を1% ギ酸の最終 V/V% に酸性化する必要が あります。4µLの10%FAを、ステップ2 の還元された溶液に添加します。静かに ボルテックスし、スピンダウンします。

 160 µL の 0.1% ギ酸を添加し、200 µL の 最終容量にします。還元されたトラスツズ マブ溶液の濃度は 250 ng/µL で、LC/MS 分析に使用できるサンプルになりました。

LC/MS 分析

LC/MS 分析は、Agilent 1290 Infinity II Bio LC システム に Agilent InfinityLab Pro iQ Plus シングル四重極 LC/MS システムを組み 合わせて実施しました (図 1)。クロマトグラ フィーによる分離のために Agilent PLRP-S カ ラム (2.1 × 50 mm, 5 μ m)を使用しました。 表 1 と表 2 に LC/MS パラメータを示します。



図 1. Agilent 1290 Infinity II Bio LC システムと Agilent Pro iQ Plus シングル四重極質量分析システム

液体クロマトグラフィー

表 1. Agilent 1290 Infinity II LC メソッド

Agilent 1290 Infinity II Bio LC システム				
カラム	Agilent PLRP-S, 2.1 × 50 mm、5 µm (部品番号 PL1912-1502)			
サンプラ温度	5℃			
移動相A	水+0.1%ギ酸			
移動相B	アセトニトリル + 0.1 % ギ酸			
流量	0.5 mL/min			
注入量	2 µL			
カラム温度	℃ C			
グラジエントプログラム (インタクト)	時間(分) %B 0 10 5 60 6 10 8 10			
グラジエントプログラム(還元)	0 5 0.1 20 8 40 8.1 70 9.1 70 9.2 5 11 5			

質量分析

表 2. MS パラメータ

Agilent Pro iQ Plus シングル四重極質量分析計			
イオン源	Agilent Jet Stream ESI イオン源		
極性	E		
タイムフィルタウィンドウ	0.1分		
ストップタイム	ポンプに従う/制限なし		
MS1 スキャン範囲	m/z 1,000 ~ 3,000 (インタクト) m/z 600 ~ 2,400 (還元)		
スキャン時間	1,500 ms		
検出器ゲイン係数	1		
フラグメンタ	275 V (インタクト) 175 V (還元)		
フラグメンタランプ?	未確認		
データ保存	プロファイル		
ガス流量	12 L/min		
ネブライザ	50 psi		
シースガス流量	11 L/min		
キャピラリー電圧	4,500 V		
ノズル電圧	2,000 V		
ガス温度	350 ℃		
シースガス温度	360 ℃		
切り替えバルブ	有効:0~1分は LC 流路を廃液ラインへ、1~8分(インタクト) または1~11分(還元)は LC 流路を MS へ		
分析後のダイバータ位置	廃液へ		

データ処理

LC/MS データは Agilent OpenLab CDS ソフ トウェア、バージョン 2.8 を使用して処理しま した。図 2 にインタクトデコンボリューション パラメータを示します。還元 mAb デコンボ リューションに対し、外部バックグラウンド時 間範囲を 3.0 ~ 4.3 分に、自動デコンボリュー ション RT ウィンドウを 4.3 ~ 5.6 分に、デコ ンボリューションされた質量範囲を 10,000 ~ 60,000 Da に、絶対ノイズスレッシュホールド を 2,000 に、相対アバンダンススレッシュホー ルドを 25 % に、MW アルゴリズムをセントロ イドに設定しました。

結果と考察

インタクトトラスツズマブの トータルイオンクロマトグラム

図 3A に 500 ng のインタクトトラスツズマブの トータルイオンクロマトグラム (TIC) を示し ます。予想どおり、TIC は約 2.6 分で溶出する 単一のピークを示しています。水性度の高い 含有物や塩が質量分析システムに混入する のを防止するために、LC/MS メソッドの最初 の1 分間、LC 流路を廃液ラインへ送ります。 1~8 分間、ダイバータバルブを切り替えて、 mAb 信号が測定される iQ Plus に直接流れる ようにしました。

インタクトトラスツズマブの未加工の MS/デコンボリューションしたスペクトル

インタクトトラスツズマブの未加工の MS/デ コンボリューションしたスペクトルをそれぞれ 図 3B と図 3C に示します。Agilent Pro iQ Plus シングル四重極質量分析のスキャン範囲 は m/z 3,000 で、インタクトトラスツズマブの多 数の電荷状態を検出し、優れたデコンボリュー ションスペクトルを得ることができました。事実、 図 3B の拡大図は、トラスツズマブの1つの 電荷状態に対する実際のピーク形状を示して います。ここで示されている相対アバンダンス は図 3C のデコンボリュートしたスペクトルで の所見と一致しており、デコンボリュートした スペクトルと、同定された5つの主なグリコ シル化ピークに対して信頼性が向上していま す。デコンボリュートしたスペクトルでゴース トピークを減少させた注目すべき処理パラ





図 2. インタクト mAb デコンボリューションに使用した、Agilent OpenLab CDS ソフトウェア、 バージョン 2.8 の処理パラメータ

メータは、OpenLab CDS MS スペクトルデー タデコンボリューションタブの基本設定に示さ れた、セット中の最小ピークのパラメータでし た (図 2)。インタクトおよび還元 mAb デコ ンボリューションに対し、この設定を 15 に設 定しました。これにより、ソフトウェアに対し、 15 以上の電荷状態を有するピークをデコン ボリューションするよう指示されます。このパ ラメータは、デコンボリュートしたスペクトル で検出されたピークの品質を向上させるため 重要です。分析対象分子の未加工の質量スペ クトルで観察された電荷状態の数に基づき、 このパラメータを調整できます。



図3. (A) Agilent Pro iQ Plus シングル四重極質量分析計で測定したインタクトトラスツズマブの TIC。(B) トラスツズマブの未加工の MS。 拡大図は 1 つの電荷状態の 拡大領域を示しています。(C) トラスツズマブのデコンボリューションした質量スペクトル、および測定質量と提示されたグリコフォーム。 表 1 および表 2 に LC/MS 設定を、 図 2 に処理パラメータを、表 3 に算出質量誤差を示します。

還元 mAb クロマトグラム、未加工の MS、デコンボリューションしたスペクトル

グリコシル化プロファイリングにさらなる側面 を追加するために、インタクト mAb をサブユ ニットへと還元し、複雑さを軽減することが可 能です。図 4A に還元トラスツズマブの TIC ク ロマトグラムを示します。軽鎖と重鎖に対する 2 つのピークを確認できます。軽鎖に対して測 定された MS スペクトル強度は重鎖の 4 倍の 高さでした。しかし、HC に対する相当な量の グリコシル化情報と、優れた MS スペクトル品 質は、Agilent Pro iQ Plus で得られました。 図 4B と図 4D の拡大図は、トラスツズマブ LC と HC の領域をそれぞれ拡大したものです。 グリコシル化部位はトラスツズマブのコンセン サス配列に一致しているため(グリコシル化 が、NXS/T モチーフに続くアスパラギン残基 で生じている)、LC は単一のピークを示すと 予想されます。アミノ酸配列によりグリカン修 飾が生じないためです。一方、トラスツズマブ HC は NXS/T モチーフを有し、そのため Fc



図 4. (A) Agilent Pro iQ Plus シングル四重極質量分析計で測定した還元トラスツズマブの TIC。トラスツズマブの軽鎖および重鎖の未加工の MS (B+D)。拡大図は、機器の 未加工のスペクトル品質を示しています。トラスツズマブの軽鎖および重鎖のデコンボリューションした質量スペクトル(C+E)。表 3 は算出質量誤差を示しています。

領域のアスパラギン残基でグリコシル化が可能 です。図 4D の拡大図は、未加工の MS スペク トルの4つのピークを示しており、これらは GO、GOF、G1F、G2F グリカンと一致します。 LC および HC のデコンボリューションしたスペ クトルをそれぞれ図 4C および図 4E に示しま す。 インタクト mAb に対する図 2 に示された パラメータと同様の処理パラメータを還元トラ スツズマブに使用しました。デコンボリューショ ンされた質量範囲を 10,000 ~ 60,000 Da に 変更し、相対アバンダンススレッシュホールドを 25%に設定したこと以外、違いはありません。 デコンボリュートしたトラスツズマブ HC のスペ クトルで注目すべきは、グリコシル化ピークの 相対強度が未加工の MS スペクトル拡大図で 観察されるものと類似しているという点です。 HC の各電荷状態はグリコシル化ピークのさま ざまな相対強度を示しますが、入力 MS スペク トルからの相対アバンダンスをデコンボリュー トしたスペクトルに関連付けすることが可能 で、それによりデコンボリューションしたスペク トルの信頼性が向上します。 還元/インタクト レベルでのグリコシル化の相対アバンダンス についてレポートできるということは重要な性 能です。これらのアバンダンスの基準値からの 偏差により、バイオ医薬品ラボの QA/QC チー ムに対してさらなる調査の実施が促されるから です。

インタクト/サブユニットレベルでの相対 アバンダンスと MS スペクトルデコンボ リューションレポート

表 4 に、インタクト/サブユニットレベルそれ ぞれにおけるトラスツズマブの測定相対アバン ダンスを示します。相対アバンダンス(%)は、 最も強い成分に対してアバンダンスを正規化 することにより計算しますが、相対定量(%) では、すべての成分の合計アバンダンスに対 する任意の成分の絶対アバンダンスを測定し ます。相対アバンダンス(%)と相対定量(%) はどちらも、製品の再現性と品質のモニタリン グに使用できます。

表 3. Agilent OpenLab CDS ソフトウェアによるデコンボリューション後の、理論上の平均分子質量と 実験で得られた分子質量の比較

Agilent Pro iQ Plus 質量分析システムでの実験で得られた測定誤差					
分子	修飾	理論上の質量(Da)	実験で得られた質量 (Da)	∆ 質量 (Da)	質量誤差 (ppm)
インタクトトラスツズマブ	G0+G0F	147,912.7	147,916.7	4.0	27
	G0F+G0F	148,058.8	148,055.1	-3.7	-25
	G0F+G1F	148,221.0	148,217.7	-3.3	-22
	G1F+G1F	148,383.1	148,374.7	-8.4	-57
	G1F+G2F	148,545.3	148,542.0	-3.3	-22
トラスツズマブ HC	GO	50,456.1	50,447.3	-8.8	-174
	G0F	50,602.2	50,597.9	-4.3	-85
	G1F	50,764.4	50,759.9	-4.5	-89
	G2F	50,926.5	50,915.4	-11.1	-218
トラスツズマブ LC	なし	23,443.3	23,440.5	-2.8	-119

表 4. (A) インタクトおよび (B) 還元レベルでのデコンボリューション後の相対アバンダンス

А

トラスツズマブインタクトの相対アバンダンス					
成分	測定質量	提示された同定	相対アバンダンス(%)	相対定量(%)	
А	148,218 Da	G0F+G1F	100	33.77	
В	148,375 Da	G1F+G1F	76.63	25.88	
С	148,055 Da	G0F+G0F	67.72	22.87	
D	148,542 Da	G1F+G2F	38.00	12.83	
E	147,917 Da	G0+G0F	13.79	4.66	

В

トラスツズマブ重鎖の相対アバンダンス				
成分	測定質量	提示された同定	相対アバンダンス(%)	相対定量(%)
А	50,598 Da	GOF	100	41.22
В	50,760 Da	G1F	94.13	38.80
С	50,915 Da	G2F	33.71	13.90
D	50,447 Da	GO	14.73	6.07

図5に、インタクトトラスツズマブの MS スペ クトルデコンボリューションソフトウェアのレ ポートを示します。レポートは OpenLab CDS ソフトウェアのレポートエディタタブで簡単に カスタマイズできます。レポートには、トレー サビリティのために、取り込みメソッドと処理 メソッド、注入量、バイアル位置、注入日に関 する詳細情報が表示されます。さらに、未加工 の MS/デコンボリューションしたスペクトルと、 検出された成分が示されます。またレポートに は、デコンボリューションされたスペクトルの成 分の相対アバンダンス(%)と相対定量(%) が一覧表示されます。

結論

LC/MS によるインタクト質量分析は、グリ コシル化の相対アバンダンスをモニタリング するために、バイオ医薬品ラボで一般的に採 用されている技術です。これらの分析は製品 の効能と安全性にとって重要です。Agilent 1290 Infinity II Bio LC システム と Agilent InfinityLab Pro iQ Plus の組み合わせは、 ルーチン分析のための堅牢なシステムを必要と するラボにとって、小型でコスト効率に優れた LC/MS 機器となります。この研究では、医薬 品のインタクトおよび還元レベルにおけるグリ コシル化ピークの相対アバンダンスをレポート するために、Pro iQ Plus LC/MS システムを 使用しました。インタクトレベルで、アジレント の高分解能精密質量分析システムで測定済み の結果と一致する相対アバンダンスにより、 5 つのグリコシル化ピークが同定されました。 トラスツズマブの重鎖に対しても相対アバンダ ンスをモニタリングし、GO、GOF、G1F、G2F グリカンと一致する 4 つのピークが同定され ました。モニタリング対象のグリコシル化の信 頼性の高い相対アバンダンスと、優れた MS スペクトルピーク形状の両方が、Agilent Pro iQ Plus シングル四重極質量分析システムの 性能を際立たせています。



図 5. インタクトトラスツズマブのMS スペクトルデコンボリューションのレポート

参考文献

- Higel, F.; Seidl, A.; Sörgel, F.; Friess, W. N-glycosylation Heterogeneity and the Influence on Structure, Function and Pharmacokinetics of Monoclonal Antibodies And Fc Fusion Proteins. Eur. J. Pharma. Biopharm. **2016**, 100, 94–100. https://doi.org/10.1016/j. ejpb.2016.01.005
- Wang, D.; Baudys, J.; Bundy, J.; Solano, M.; Keppel, T.; Barr, J. Comprehensive Analysis of the Glycan Component of SARS-CoV-2 Spike Proteins Using Signature Ions-Triggered Electron-Transfer/Higher-Energy Collisional Dissociation (EThcD) Mass spectrometry. Anal. Chem. **2020**, 92, 21, 14730-14739. https://doi.org/10.1021/acs. analchem.0c03301

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。

DE-006222

アジレント・テクノロジー株式会社 © Agilent Technologies, Inc. 2025 Printed in Japan, May 13, 2025 5994-8345JAJP

