バイオ医薬品

# 複数の GLP-1 アナログの包括的な特性解析

Agilent 6545XT AdvanceBio LC/Q-TOF と電子捕獲解離 および ExDViewer ソフトウェアの使用

## 著者

Rachel Franklin, Mike Hare, Thomas Walker, and Joseph Meeuwsen Agilent Technologies, Inc.

## 概要

GLP-1 アナログは、グルカゴン様ペプチド-1 (GLP-1)と呼ばれる天然由来のペプチドホル モンに似た合成ペプチドです。GLP-1 アナログは糖尿病と肥満を管理できる可能性がある ため、その開発は重要です。アジレントの高分解能 LC/Q-TOF を使用して合成ペプチドを 特性解析すると、GLP-1 アナログに関する重要な品質情報を得ることができます。LC/MS ワークフローに電子ベースのフラグメンテーションを組み込むことは、タンパク質とペプ チドの配列と修飾を分析するための効果的な手法です。このアプリケーションノートでは、 Agilent 6545XT AdvanceBio LC/Q-TOF システムと電子捕獲解離用の Agilent ExD セル による 3 種類の GLP-1 アナログ(リラグルチド、セマグルチド、チルゼパチド)の特性解析 について説明します。Agilent ExDViewer は、タンパク質やペプチドから得られるあらゆる 種類のフラグメンテーションデータの解析に適した使いやすいソフトウェアツールです。この アプリケーションノートでは、ExDViewer のターゲットデコンボリューションワークフローを 使用して詳細な配列分析とカスタム修飾の特性解析を実行し、強力なツールでフラグメン テーションの傾向とデータ品質を視覚化します。

#### 概要

GLP-1 は GLP-1 受容体と結合するインスリン刺激ホルモンで す。<sup>1、2</sup> 2019 年に最初の GLP-1 アナログが II 型糖尿病向け に米国食品医薬品局 (FDA) によって承認され、2020 年には GLP-1 アナログが肥満管理向けに承認されました。<sup>3、4</sup> GLP-1 アナログの使用は増えているため、高精度な質量分析 (MS) メ ソッドを構造特性解析、純度、偽造品の分析に使用することが重 要です。

電子捕獲解離(ECD)による合成ペプチドのLC/MS/MS分析により、配列、修飾、純度の詳細情報を得ることができます。 ECDは衝突誘起解離(CID)とは異なり、GLP-1アナログ中の 脂肪酸修飾などの修飾を保持する「緩やかな」フラグメンテー ションアプローチと見なされています。このため、ペプチドバック ボーンを効果的に切断しながら、マッピングや特性解析を実行で きます(図1)。また、ECDによるアミノ酸側鎖の二次的フラグメ ンテーションにより、配列割り当てのさらなる証拠を入手し、アス パラギン酸やイソアスパラギン酸などのアミノ酸異性体を調査で きます。<sup>5.6</sup> Agilent ExD セルと Agilent ExDViewer ソフトウェア は、複雑な構造を持つ合成ペプチドの詳細な特性解析に最適で す。

このアプリケーションノートでは、3 種類の GLP-1 アナログ(セ マグルチド、リラグルチド、チルゼパチド)の LC/MS/MS 分析に ついて説明します。この分析は、6545XT LC/Q-TOF と ExD セ ルを使用し、ECD を有効にして実行しました。ExDViewer を使 用して、ターゲットデコンボリューションワークフローを用いた MS/MS フラグメンテーションスペクトル分析を実施しました。こ のワークフローでは、電子フラグメンテーション独自の特性(水 素移動や側鎖フラグメンテーションなど)が考慮されます。<sup>7</sup>また ExDViewer には、ECD と CID の両方の実験でフラグメンテー ションの傾向を視覚化し、データ品質を把握するための便利な ツールも含まれています。これらのメソッドでは、ExD のハード ウェアとソフトウェアを組み合わせることで、非標準の残留物を 含むペプチド配列の高速で詳細な特性解析と、カスタム修飾の 特性解析を実行できました。



図 1.アミノ酸位置 2 に非標準アミノ酸である 2-アミノイソ酪酸(文字 B で表示)が 含まれるセマグルチドの構造。リジン 20 は、ビス-アミノジエトキシセチル(2xAdo) とガンマグルタミン酸リンカーにより結合している C18 脂肪酸によって修飾されて います。リラグルチドとチルゼパチドのリジン修飾は構造が似ており、リンカーと脂 肪酸基が異なります。

## 実験方法

### 試薬と標準試料

- Agilent ESI-L チューニング混合液(部品番号 G1969-85000)、2.5 µg/mL のメリチン(部品番号 G1997-85001) を含む
- ギ酸、99.0%以上、Optima LC/MS グレード、カタログ番号
  A-117-50、Fisher Chemical、(部品番号 A-117-50)
- アセトニトリル、99.9 % 以上、LC/MS グレード、Supelco OmniSolv (部品番号 AX0156-6)
- リラグルチド、Sigma (部品番号 SML3925)
- セマグルチド、AstaTech (部品番号 AT35750)
- チルゼパチド、AstaTech (部品番号 AT40456)

#### サンプル前処理

各サンプルは、0.1 % のギ酸を含む 15 % のアセトニトリル 10 μM で前処理しました。 LC 実験用に、等量の 3 種類のサンプ ルを混合しました。

#### HPLC カラム

Agilent AdvanceBio ペプチドマッピング 120 Å、2.1 × 150 mm、2.7 µm (部品番号 653750-902)

## 装置構成

- Agilent 1290 Infinity II Bio LC システム
  - Agilent 1290 Infinity II ハイスピードポンプ (G7120A)
  - Agilent 1290 Infinity II マルチサンプラ (G7167B)
  - Agilent 1290 Infinity II カラムコンパートメント (G7116B)
- Agilent 6545XT AdvanceBio LC/Q-TOF MS と Agilent ExD セル

## ソフトウェア

- Agilent MassHunter Acquisition ソフトウェア(LC/Q-TOF 用)、バージョン 11.0
- Agilent ExDControl ソフトウェア、バージョン 3.6
- Agilent ExDViewer ソフトウェア、バージョン 4.6.12

#### 液体クロマトグラフィーメソッド

このアプリケーションノートに記載されている LC メソッドを、ほ とんどの ECD 実験に使用しました。ただし、ExD セルのチュー ニングと、荷電状態に依存するフラグメンテーションの調査には、 直接注入を使用しました。直接注入は、500 µL シリンジを使用 して、注入速度 10 ~ 20 µL/min で実行しました。PEEK チュー ブを使用して、注入シリンジから Agilent Dual AJS ソースのネブ ライザ注入口にサンプルを移しました。図 2 に、3 種類の GLP-1 アナログすべての混合物のクロマトグラフィーによる分離を示し ます。

Agilent 1290 Infinity II Bio LC システム				
カラム	Agilent AdvanceBio ペプチドマッピング 120 Å、 2.1 × 150 mm、2.7 µm			
移動相A	LC/MS グレードの水 + 0.1 % ギ酸			
移動相 B	アセトニトリル + 0.1% ギ酸			
流量	0.400 mL/min			
注入量	10 µL			
カラム温度	60 °C			
グラジエントプログラム	時間(分) 0	%B 20		
	20	60		
	20.10	80		
	21.90	80		
	22	20		

表 1. GLP-1 アナログに使用した液体クロマトグラフィーメソッド



図 2. GLP-1 アナログの混合物のクロマトグラム。ピーク ID は次のとおりです。 (A) セマグルチド、(B) リラグルチド、(C) チルゼパチド

#### 質量分析メソッド

#### 表 2. LC/Q-TOF パラメータ

Agilent 6545XT Q-TOF MS システム		
イオン源	Agilent Dual Jet Stream ESI イオン源	
極性	Æ	
ガス温度	325 °C	
乾燥ガス流量	10 L/min	
ネブライザ	35 psi	
シースガス温度	275 °C	
シースガス流量	12 L/min	
キャピラリ電圧	3,500 V	
ノズル電圧	2,000 V	
フラグメンタ	175 V	
スキマ電圧	75 V	
取り込みレート	2 スペクトル/秒	
取り込みモード	拡張ダイナミックレンジ(2 GHz)	
四重極選択幅	中 (4 amu)	
MS スペクトル範囲	m/z 120 ~ 3,200	
MS/MS スペクトル範囲	m/z 120 ~ 3,200	

#### 表3. 自動 MS/MS データ取り込みパラメータ

パラメータ	設定値
	4+ プリカーサ用の推奨リスト:
プリカーサの選択	- m/z 938.2 (リラグルチド)
	- m/z 1,028.7(セマグルチド)
	- m/z 1,203.6 (チルゼパチド)
強度スレッシュホールド(Abs)	2,000
強度スレッシュホールド(Rel)	0.01 %
質量誤差許容範囲	20 ppm
静的排除	m/z 100 ~ 400
同位体モデル	ペプチド
推奨イオンリストのみを使用	無効

## ターゲット ECD MS ワークフロー

ExD セルレンズと内部フィラメントは、ExDControl ソフトウェアバージョ ン 3.6 で制御しました。ExDControl は MassHunter Acquisition ソフ トウェアと共に、別個のアプリケーションとして動作します。最適なフィラ メント電流 (2.55 A) をフィラメントのオートチューンで決定してから、手 動で調整してメリチンの ECD フラグメントを最適化しました。インタクト イオンの MS1 透過率用のレンズプロファイルは、チューンミックスイオン でオートチューンして決定しました。メリチンの ECD フラグメントイオン ピークを使用して、MS2 の ECD に使用するレンズプロファイルをオート チューンしました。LC 実験のために、表 2 と表 3 のパラメータを使用し て、MassHunter Acquisition ソフトウェアで自動 MS/MS 取り込みメ ソッドを設定しました。直接注入実験のために、ターゲット MS/MS 取り 込みを使用して、個々の荷電状態を調査しました。

#### データ解析

GLP-1 アナログの分子量は、ExDViewer の MS1 デコンボリューション ワークフローを使用して測定しました。デコンボリューションのために未 加工の.d ファイルを ExDViewer に直接読み込み、デフォルトパラメータ を使用しました。分離ピークごとに MS1 スペクトルを平均化しました。 スペクトルウィンドウのスペクトルビューを切り替えて、放電スペクトルを 生成しました。ファイルは.svg ファイルとして保存し、図要素のカスタマ イズのために Microsoft PowerPoint にインポートしました。

配列と修飾の分析には、ExDViewer のターゲットデコンボリューション を使用しました。まず、ターゲットエディタで GLP-1 アナログの配列を定 義し、MS/MS 分析結果に対して検索しました。 フラグメント照合には デフォルトのパラメータを使用しました。ただし、リジン修飾の特性解析 は除きます。修飾分析では、イオンスコアのスレッシュホールドを 10 に 上げて、フラグメント照合を含む最高品質の修飾のみを観察しました。



図 3. GLP-1 アナログ合成ペプチドの LC/MS/MS 分析のワークフロー

# 結果と考察

## ExDViewer による同位体分離されたスペクトルデータの デコンボリューション

Agilent LC/MS/MS ワークフローを用いた GLP-1 アナログの分析に より、1 回の実験で分子量、信号の不均一性、配列を測定できます。 ExDViewer のノンターゲットデコンボリューションアルゴリズムが、(通常 はペプチド分析中に得られる) 同位体分離されたスペクトルからの信号 を効果的に処理します。MS1 デコンボリューションにより、合成ペプチド の分子量を測定できます。 図 4 に、図 2 の混合物からの各 GLP-1 アナログの MS1 とデコンボ リューションしたスペクトルを示します。デコンボリューションしたスペク トルによって、GLP-1 アナログに対応する各分離ピークからの一次信号 と、低アバンダンスの塩が付加されたいくつかのピークが現れます(Na<sup>+</sup>/ K<sup>+</sup>)。ExDViewer による同位体分離された信号のデコンボリューション は、Agilent MassHunter BioConfirm のより大きいタンパク質イオンの 最大エントロピーデコンボリューションを補完するものです。



図 4.LC 分離した GLP-1 アナログの MS1 スペクトル (左)と対応するデコンボリュートしたスペクトル (右): (A) セマグルチド、(B) リラグルチド、(C) チルゼパチド。 モノアイソトピック質量はデコンボリュートした質量スペクトルでアノテーション済みです。

#### 非標準アミノ酸による配列分析

ECD によるガス相のフラグメンテーションによりペプチドバックボーンが 正確に切断され、ExDViewer で容易に解釈できる明確で再現性の高い フラグメンテーションスペクトルが生成されます。セマグルチドとチルゼパ チドには非標準アミノ酸が含まれています。これらは ExDViewer 内の構 成要素 (Building Block) エディタでカスタムアミノ酸と定義される場合 があります。今回の分析では、非標準アミノ酸を 2-アミノイソ酪酸と定義 し、省略コードとして B の 1 文字を使用します。2 Hz で収集したシング ルターゲット MS/MS スペクトルを使用して、リラグルチド、セマグルチド、 チルゼパチドの 4+ プリカーサの 100 % の配列カバーを取得しました。 CID と ECD の複数のフラグメントイオンタイプ (a、b、c、d と w、y、z) が検出されました。これが配列確認の補完的な証拠となります (図 5)。 ExD セルを使用すると、一般的に豊富なアミノ酸側鎖フラグメントが生成



図 5. 配列カバーマップ:(A)セマグルチド、(B)リラグルチド、(C)チルゼパチド。非標準の B アミノ酸は黄色で強調表示しており、修飾済みリジンは赤文字で示しています。 c/z タイプと b/y タイプの相補的イオンがいくつか同定されました。Agilent ExDViewer フラグメント照合にはデフォルトのパラメータを使用しました。

され、ロイシンやイソロイシンなどの異性体アミノ酸に関する詳細情報を 取得できます。ECD ラジカルイオンからのアミノ酸側鎖の二次的フラグメ ンテーションは、W イオンと呼ばれます。<sup>4</sup> ExDViewer は側鎖の証拠を 自動的にアノテーションするため、アミノ酸異性体の分析が容易になりま す。<sup>7</sup> イソアスパラギン酸形成に固有のイオンはタンパク質品質の評価用 の重要な修飾であり、これもアノテーションされます。<sup>6.7</sup>図6では W<sub>23</sub><sup>2+</sup> イオンがアノテーションされており、この位置のアスパラギン酸のさらなる 証拠となります。W<sub>23</sub><sup>2+</sup> イオンにカーソルを合わせると情報ツールチップ が開き、水素移動率、質量、側鎖喪失の式、スコア、一致する同位体の 数が表示されます。



図 6.リラグルチドの代表的なフラグメンテーションスペクトル。アスパラギン酸やイソアスパラギン酸などの異性体アミノ酸の自動的にアノテーションされたフラグメントの証拠を示しています。アスパラギン酸は、Z23<sup>2+</sup> イオンからの CHO2 基の喪失により特定されます。

#### ECD と CID によるフラグメンテーションの傾向の視覚化

ExDViewer で自動生成されるプロットに、さまざまな実験条件でのフラ グメンテーションの傾向が表示されます。フラグメンテーションの傾向を 視覚化すると、データ品質を評価できます。例えばイオン強度のグラフに は、各種イオン(内部イオンや割り当てられないイオンなど)で検出され る信号が表示されます。これらのグラフを使用して、コリジョンエネルギー などのメソッドパラメータの適合性を評価することもできます。図7 では イオン強度グラフを使用して、ECD または CID と ECD の組み合わせを 用いた場合の、荷電状態に依存するセマグルチドのフラグメンテーション パターンを示しています。 コリジョンエネルギーがゼロの場合、5+ および 4+ セマグルチドプリカー サの 100 % 配列確認を ECD で問題なく実行できました。ただし 3+ プリ カーサの場合、ECD だけではグラフメンテーションが困難でした。35 V のコリジョンエネルギーを追加すると、3+ プリカーサの配列カバーが 57 % から 93 % に改善されました (図 7C と図 7F)。注目すべきなのは、 3+ セマグルチドのフラグメントイオンアバンダンスの合計が、CID の有無 にかかわらず同等であることです。つまり、CID 追加時の配列カバーの増 大は、フラグメントイオンアバンダンス全般の増加ではなく、別のフラグメ ント経路を反映しています。したがって、ECD と低レベルの活性化を使用



図 7. セマグルチド 3+、4+、5+ プリカーサイオンのフラグメント経路。A、B、C のパネルは、ECD のみの条件下でのイオン強度を示しています。D、E、F のパネルは、ECD と 35 V の衝突活性化の組み合わせ (ECD+CID) によるイオン強度を示しています。これらのデータセットの分析において、内部フラグメントイオンは考慮していません。

すれば、低電荷ペプチドでもより完全な配列カバーを達成できます。ただし、過大なコリジョンエネルギーを加えると、割り当てられないイオン信号の割合が高くなります。これはおそらく過剰なフラグメンテーションによるものです(図7D)。

## GLP-1 アナログ修飾の特性解析

GLP-1 アナログの新しい派生物が作られているため、その修飾の特性解析に適したメソッドを用意することが重要です。今回はフラグメント分析 に ExDViewer を使用して、3 種類の GLP-1 アナログの合成修飾の部 位と構造を特定しました。GLP-1 アナログ修飾のカスタムケミストリは、 ExDViewer の修飾エディタ (Modifications Editor) で定義しました。 ExDViewer のフラグメント照合アルゴリズムは、*m/z* と強度一致、およ び周囲のノイズレベルに基づいて各同位体クラスタとイオンスコアを割り 当てます。デフォルトのイオンスコア設定は、多くの配列分析アプリケー ションに使用できます。ただし、修飾の特性解析には、修飾位置を特定す るための厳密なスペクトル品質が必要です。図8に、チルゼパチドで検 出された、修飾を含むいくつかの ECDイオンの例を示します。この図から、 フラグメンテーションスペクトルで見られるイオン照合品質の範囲がわか ります。



図8.さまざまなイオンスコアのチルゼパチドの代表的なフラグメント。各イオンには、m/z と予想される強度一致、および周囲のノイズレベルに基づいてスコアが割り当てられます。(A) チルゼパチドの配列カバーマップ。スコアが5以上のイオンのみを考慮しています。(B ~ E) フラグメントとその対応するイオンスコアを含む修飾の例。修飾済みのリジンは赤文字で示しています。C20<sup>24</sup> イオンは部位固有の局在化フラグメントです。その他の強調表示されているイオンは、ペプチドバックボーンのさまざまな位置でフラグメント 化された修飾を含むペプチドです。

ECD は衝突解離とは異なり、重要かつ不安定な化学修飾を保持しなが らペプチドバックボーンをフラグメント化する、緩やかな手法です。ECD や CID で分析した GLP-1 アナログのイオンスコアを比較すると、ECD フ ラグメンテーションでは CID よりも高品質な、修飾を含むフラグメントが 生成されることがわかります(図 9)。修飾の両側の高品質な照合によ り、局在診断の信頼性が向上します。多くの場合、ECD フラグメンテーショ ンだけで、コリジョンエネルギーを最適化しなくても詳細な構造情報を取 得できるため、配列分析の作業時間を短縮できます。



**図 9.** ECD または CID の実験で同定された、イオンスコアが 10 以上のチルゼパチドフラグメント。ECD では、特に修飾の N 末端側で、より高品質な修飾を含むフラグメントを 得ることができました。ECD では、相補的イオンタイプにより、イオン化列がより完全なものになりました。この例外はプロリンが豊富な領域です。ECD はプロリンの N 末端側 を切断しないためです。

# 結論

このアプリケーションノートでは、3 種類の合成 GLP-1 アナログでの電子 解離による包括的な配列分析と局在診断について説明しています。フラ グメント分析には Agilent 6545XT AdvanceBio LC/Q-TOF と Agilent ExDViewer を使用しました。アミノ酸側鎖のフラグメンテーションにより、 アスパラギン酸やイソアスパラギン酸などのアミノ酸異性体を同定するた めの、さらなる証拠を得ることができます。ECD データからの高品質なフ ラグメント照合により、修飾の特性解析と局在診断を効率化できます。こ れはコリジョンエネルギーの最適化が不要になるためです。ExDViewer は、ECD と CID の実験でフラグメント経路とデータ品質を視覚化するの に最適なツールです。このアプリケーションノートで紹介したメソッドは、 GLP-1 アナログの分析の枠組みとなるものです。また、ECD を使用して も使用してください。

# 参考文献

- Andersen, A.; Lund, A.; Knop, F. K.; Vilsbøll, T. Glucagon-Like Peptide 1 in Health and Disease. *Nat. Rev. Endocrinol.* **2018**, 14(7), 390-403.
- Holst, J. J. The Physiology of Glucagon-Like Peptide 1. *Physiol. Rev.* 2007, 87(4), 1409–1439.
- 3. U.S. Food and Drug Administration FDA News Release. FDA Approves First Oral GLP-1 Treatment for Type 2 Diabetes. https://www.fda.gov/news-events/press-announcements/ fda-approves-first-oral-glp-1-treatment-type-2-diabetes (accessed 2019).
- 4. U.S. Food and Drug Administration FDA News Release. FDA Approves Weight Management Drug for Patients Aged 12 and Older. https://www.fda.gov/drugs/news-events-humandrugs/fda-approves-treatment-adults-alzheimers-disease (accessed 2020).
- Beckman, J. S.; Voinov, V. G.; Hare, M.; Sturgeon, D.; Vasil'ev, Y.; Oppenheimer, D.; Shaw, J. B.; Wu, S.; Glaskin, R.; Klein, C.; et al. Improved Protein and PTM Characterization with a Practical Electron-Based Fragmentation on Q-TOF Instruments. J. Am. Soc.Mass Spectrom. 2021, 32(8), 2081-2091.
- Wang, J.; Mukherjee, S.; Zubarev, R. A. Isoaspartate and Neurodegeneration. *Aging* (Albany NY), **2022**, 14(22), 8882.
- Identification of Amino Acid Isomers Using Electron Capture Dissociation in the Agilent 6545XT AdvanceBio LC/Q-TOF System. Agilent Technologies application note, publication number 5994-7506EN, 2024.

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。

#### RA250218.98

アジレント・テクノロジー株式会社 © Agilent Technologies, Inc. 2025 Printed in Japan, April 14, 2025 5994-7994JAJP

