

# 高分離能サンプリング 2D-LC による 医薬品不純物分析

API ピークに隠れた低濃度不純物の検出

# アプリケーションノート

低分子医薬品およびジェネリック医薬品

# 著者

Susanne Stephan and Sonja Krieger Agilent Technologies, Inc. Waldbronn, Germany

# 概要

医薬品有効成分 (API) に対する低濃度の不純物の分析は原薬の品質管理において重要です。不純物 が API と構造的に類似し、濃度差が大きい場合、クロマトグラフィーによる分離や検出が困難な場合 があります。

Agilent 1290 Infinity II 2D-LC ソリューションでは、コンプリヘンシブ (LCxLC) とマルチハートカット (MHC) と高分離能サンプリング 2D-LC (HiRes 2D-LC) との間の切り替えが容易に行えます。このアプリケー ションでは、近接して溶出する 2 つの化合物を HiRes 2D-LC によって分離します。このうちの 1 つは、 きわめて低い濃度で存在し、もう一方の高濃度の化合物に隠れています。クロロジフルオロ安息香 酸および脱アミドインスリンをそれぞれ標準物質および実際のサンプルで分析しました。





# はじめに

原薬の純度分析では、品質および患者の安 全がきわめて重要です<sup>1</sup>。ICH ガイドラインの 03A(R2) によると、新しい原薬内の不純物は API に対して、0.05 %<sup>2</sup> のスレッシュホールドを 超えると報告が必要になります。この濃度の 不純物を分離して検出するのは困難な場合 があります。特に、不純物が医薬品有効成分 (API) と構造的に類似している場合です。

このアプリケーションノートでは、不純物の分 離と検出に高分離能 (HiRes) サンプリング 2D-LC を用いて、相対濃度 100 % と 0.05 % で存 在するクロロジフルオロ安息香酸の各種異 性体の分析について説明します。<sup>2</sup>D 検出器と して、Agilent 1290 Infinity II ハイダイナミックレ ンジダイオードアレイ検出器 (HDR-DAD) 不純 物分析ソリューションを用いたこのアプリケー ションで、精度と真度をさらに向上することが できました。脱アミドインスリンは実際のサン プルとして、HiRes サンプリング 2D-LC/MS を 用いて分析し、構造的に類似した不純物の検 出を示します。2D-LC のさらなる利点として、 <sup>2</sup>D 分離が脱塩ステップとなり、MS イオン源に おいて<sup>1</sup>D 緩衝液からの高濃度の塩によって 生じるイオン抑制を回避することができます3。 Agilent MassHunter O<sup>2</sup>D Chromatogram Creator を用いて MassHunter の 2D-LC/MS データの表 示と解析が可能です。

図1に示したように、HiRes サンプリング2D-LC では、選択した時間範囲にわたっていくつ かの小さいフラクションを収集することでター ゲット化合物を測定でき、<sup>1</sup>D クロマトグラムの 1 つのピーク領域全体が対象となるように設 定できます。各カットがサンプリングループ内 で貯められ、すべてのカットが連続して二次元 目で分析されます。このモードで、選択した化 合物の総量が二次元目に確実に送られ分析 されます。このように、選択した共溶出化合物 を二次元目での分析用として、HiRes サンプリ ングプロセスに導入できます。こうすることに より、技術概要<sup>4</sup>に従って、信頼性の高い定量 が可能になりました。

# 実験方法

#### 装置

Agilent 1290 Infinity II 2D-LC ソリューションは、 次のモジュールで構成しました。

- ・ 2 台の 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)、Infinity II サンプル冷却システム (オプション #100) を搭載
- 2 台の Agilent 1290 Infinity II マルチカラム サーモスタット (G7116B)
- 3 台の Agilent 1290 Infinity II ダイオード アレイ検出器 (G7117B)、光路長 3.7 mm、 10 mm、60 mm の Max-Light カートリッジ セル (G4212-60008) を搭載
- Agilent 1290 Infinity バルブドライブ (G1170A)、2 ポジション/4 ポートデュオ バルブ (2D-LC バルブヘッド、G4236A) を 搭載
- 2 台の Agilent 1290 Infinity バルブドライブ (G1170A)、40 µL ループ付きマルチハート カットバルブ (G4242-64000) を搭載

脱アミドインスリンは、Agilent 1290 Infinity II 2D-LC ソリューションとデュアル Agilent Jet Stream ESI ソースを備えた Agilent 6545 0-TOF を組み 合わせて分析しました。以前のアプリケーショ ンノート<sup>5</sup>で推奨したように、<sup>1</sup>D 緩衝液中の大 量の塩がイオン源に入るのを避けるために、 タイムテーブルを用いて MS ダイバータバルブ を切り替え、<sup>2</sup>D 分析の最初の 1 分間は廃液へ 流れるようにします。

#### ソフトウェア

- Agilent OpenLab CDS ChemStation Edition Rev. C.01.07 SR2 [255]、Agilent 2D-LC ソフトウェア (バージョン A.01.03 [025])、 および Agilent HDR-DAD ChemStation アドオン (製品バージョンA.01.01 [015]) 付属
- Agilent MassHunter ワークステーション ソフトウェア LC/MS データ取り込み、 バージョン B.08.00、ビルド 8.00.8026.0
- Agilent MassHunter ワークステーションソ フトウェア Qualitative Analysis、バージョン B.07.00、ビルド 7.0.7024.0
- ・ MassHunter Rev. 1.0.15 の Agilent <sup>2</sup>D Chromatogram Creator



図 1. 高分離能サンプリング 2D-LC の説明図

## 試薬

すべての溶媒は LC グレードを使用しました。 アセトニトリルおよび硫酸ナトリウムは、Merck (ダルムシュタット、ドイツ)から購入しました。 ギ酸、リン酸、二塩基性リン酸ナトリウムと ー塩基性リン酸アンモニウムは Sigma-Aldrich (シュタインハイム、ドイツ)から入手しました。 超純水は、0.22 µm メンブレンユースポイント カートリッジ (Millipak、EMD Millipore、ビレリカ、 マサチューセッツ州、米国)を装備した Milli-O Integral システムで精製しました。標準物質の 5-クロロ-2,4-ジフルオロ安息香酸、3-クロロ -2,4-ジフルオロ安息香酸、2-クロロ-4,5-ジフル オロ安息香酸、ウシ膵臓由来のインスリンは Sigma-Aldrich (シュタインハイム、ドイツ)から 購入しました。

#### サンプル

クロロジフルオロ安息香酸の原液は、 水/アセトニトリル (50/50)を用いて1 mg/mLの 濃度に調製しました。異なる濃度比の3つの 化合物の混合物を原液から調製しました。

1 mg/mL のインスリンは、25 mM リン酸ナトリ ウム緩衝液 (pH 2.55) に溶解しました。脱アミ ド化のために、インスリン溶液を pH 9 に調整 し、室温で12 時間静置しました。

## クロロジフルオロ安息香酸の分析に使用した高分離能サンプリング 2D-LC メソッド

	2D					
	DAD. 光路長 10 mm の Max-Light	<sup>2</sup> D				
	カートリッジャル付き	HDR-DAD ソリューション				
カラム						
一次元日	Agilent ZORBAX Eclinse Plus C18 RBHD					
八九日	$21 \times 100 \text{ mm}, 1.8 \text{ µm} (\text{p/n} 959758-902)$					
- 次元日	Agilent ZOBBAX Eclinse Plus PAH					
_,0001	$2.1 \times 100 \text{ mm}$ 1.8 $\mu$ m (p/n 959764-918)					
<sup>1</sup> D ポンプ						
溶媒A	水 + 0.1 % リン酸					
	アセトニトリル					
流量	0.2 mL/min					
グラジエント	0 分 – 30 %B					
	10 分-30 %B					
	12 分-80 %B					
	14 分-80 %B					
	15 分-30 %B					
<sup>2</sup> D ポンプ						
溶媒A	水 + 0.1 % リン酸					
溶媒 B	アセトニトリル					
流量	0.5 mL/min					
グラジエント	0 分-20 %B					
	4 分-25 %B					
	5 分-25 %B					
<sup>2</sup> D グラジエントストップタイム	5.00 分					
<sup>2</sup> D サイクル時間	6.00 分					
ストップタイム	15 分					
高分離能サンプリング						
タイムベース	6.54 分	7.83 分				
サンプリング時間	7秒	9秒				
カット数	9	10				
マルチカラムサーモスタット						
一次元目	30 ° C					
二次元目	30 ° C					
マルチサンプラ						
注入量	1 µL	5 µL				
ニードル洗浄	メタノール/水 (50/50) で 10 秒					
<sup>1</sup> D ダイオードアレイ検出器						
波長	210 nm/4 nm、リファレンス 395 nm/10 nm					
データレート	40 Hz					
<sup>2</sup> D 検出器						
	DAD (10 mm セル)	HDR-DAD				
波長	210 nm/4 nm、リファレンス 395 nm/10 nm					
データレート	40 Hz					

# 脱アミドインスリンの分析に使用した高分離能サンプリング 2D-LC メソッド (変更)<sup>3</sup>

パラメータ	設定値			
カラム				
一次元目	Agilent Poroshell 120 EC-C18			
	$2.1 \times 150 \text{ mm}_{\circ}$ 2.7 $\mu\text{m}$ (p/n 693775-902)			
二次元目	Agilent ZORBAX Bonus RP、			
	$2.1 \times 50 \text{ mm}$ 1.8 $\mu\text{m}$ (p/n 857768-901)			
<sup>1</sup> D ポンプ				
溶媒A	$60 \text{ mM } \text{Na}_2\text{SO}_4 + 40 \text{ mM } \text{NH}_4\text{H}_2\text{PO}_{4\text{\tiny $n$}} \text{ pH } 2.21$			
溶媒 B	アセトニトリル/水 (80/20)			
流量	0.3 mL/min			
グラジエント	0分-29%B			
	28 分 - 43 %B			
	29 分 - 90 %B			
	30 分 - 90 %B			
	31 分 - 29 %B			
<sup>2</sup> D ポンプ				
溶媒 A	水 + 0.1 % ギ酸			
溶媒 B	アセトニトリル/水 (80/20) + 0.1 % ギ酸			
流量	0.4 mL/min			
グラジエント	0分-6.25%B			
	0.1 分 - 23% B			
	2.6 分 - 35% B			
	2.7 分 - 90% B			
	2.8 分 - 90% B			
	2.9 分 - 6.25% B			
<sup>2</sup> D グラジエントストップタイム	3.00 分			
<sup>2</sup> D サイクル時間	5.00 分			
ストップタイム	38 分			
高分離能サンプリング				
タイムベース	15.64 分			
サンプリング時間	4秒			

パラメータ	設定値
カット数	10
マルチカラムサーモスタット	
一次元目	40 ° C
二次元目	40 ° C
マルチサンプラ	
注入量	1 μL
ニードル洗浄	メタノール/水 (50/50) で 10 秒
<sup>1</sup> D ダイオードアレイ検出器	
波長	195 nm/4 nm、リファレンス 395 nm/100 nm
データレート	40 Hz
<sup>2</sup> D ダイオードアレイ検出器	
波長	195 nm/4 nm、リファレンス 395 nm/100 nm
データレート	40 Hz
MS パラメータ	
モード	ポジティブ
ガス温度	200 ° C
ガス流量	13 L/min
ネブライザ	35 psig
シースガス温度	375 ° C
シースガス流量	12 L/min
VCap	2,500 V
ノズル電圧	300 V
フラグメンタ	175 V
スキマ電圧	65 V
Oct 1 RF Vpp	750 V
質量範囲	100 ~ 3,200 m/z
採取レート	2スペクトル/秒

# MS ダイバータバルブのタイムテーブル

時間セグメント 番号	開始時間 (分)	ダイバータバルブ 位置
1	0	MS
2	16.02	廃液
3	17.02	MS
4	21.02	廃液
5	22.02	MS
6	26.02	廃液
7	27.02	MS
8	31.02	廃液
9	32.02	MS
10	36.02	廃液
11	37.02	MS
12	41.02	廃液

時間セグメント 番号	開始時間 (分)	ダイバータバルブ 位置
13	42.02	MS
14	46.02	廃液
15	47.02	MS
16	51.02	廃液
17	52.02	MS
18	56.02	廃液
19	57.02	MS
20	61.02	廃液
21	62.02	MS
22	66.02	廃液
23	67.02	MS

## 高分離能サンプリング 2D-LC のための メソッド設定

Agilent 1290 Infinity II 2D-LC ソリューションによ る高分離能サンプリング 2D-LC を実施しまし た。バルブ構成 (図 2) は 12 のサンプリングルー プを保持する2個のマルチハートカットバルブ に接続された 2 ポジション/4 ポートデュオバ ルブから成ります。このセットアップにより、最 大10の連続するカットをサンプリングして分 析するまで貯めておくことができます。HiRes サ ンプリング 2D-LC では、サンプル損失を防ぐた めに80%以下のループ充填を推奨していま す。図3に、クロロジフルオロ安息香酸の分 析において<sup>2</sup>D ポンプで使用するメソッド設定 を示します。最初に、サンプルの 1D-LC 分離を 実行し、クロマトグラムをプレビューウィンド ウ内にリファレンス信号としてロードしました。 ターゲットピークについて、クロロジフルオロ 安息香酸のピーク幅全体をカバーする9つの カットによって時間ベースで HiRes サンプリン グを設定しました。提供された<sup>1</sup>D条件では、 7 秒のサンプリング時間は 58% のループ充 填に相当します。<sup>2</sup>D 検出器として Agilent 1290 Infinity II HDR-DAD 不純物分析ソリューションを 用いて注入量を増やすと<sup>1</sup>D ピーク幅全体をカ バーするには 10 カット必要となり、サンプリン グ時間が9秒と長くなります。これは75%の ループ充填に相当します。脱アミドインスリン の HiRes サンプリングでも同じ手順でメソッド を設定します。10 カットはサンプリング時間 4 秒で、50%のループ充填に相当します。



図 2.12 のサンプリングループを保持する Agilent 1290 Infinity II 2D-LC のバルブ構成



図 3. <sup>2</sup>D ポンプのメソッド設定

## 結果と考察

#### クロロジフルオロ安息香酸の分析

異なる濃度比の3 つの異性体、2-クロロ -4,5-ジフルオロ安息香酸(化合物1)、5-クロ ロ-2,4-ジフルオロ安息香酸(化合物2)、3-クロ ロ-2,4-ジフルオロ安息香酸(化合物3、構造は 図4)をHiResサンプリング2D-LC分析におい て、<sup>2</sup>D検出器として光路長10mmのMax-Light カートリッジセル付きAgilent 1290 Infinity II DAD を用いて分析しました。



図 4. 分析した 3 つの クロロジフルオロ安息香酸の化学構造

図 5 に、2 つの混合物の<sup>1</sup>D クロマトグラムを 示します。3 つすべての化合物を同じ低濃度で 含む混合物(青)と、化合物3を相対濃度100 %、化合物1と2をそれぞれ相対濃度 0.05% で含む混合物(赤)です。青のクロマトグラム では3つのピークが分離しており、赤いクロ マトグラムでは化合物 2 (相対濃度 0.05 %) は 化合物3(相対濃度100%)の下に隠れ、分離 ピークとして検出されませんでした。これらの 化合物の 2D-LC 分析で、HiRes サンプリングを 用いてピーク3の幅全体をサンプリングし、続 いて<sup>2</sup>D カラムで 9 つのカットを連続して分析 しました。図 6 に、<sup>1</sup>D ピークのサンプリングス キームと続いて得られたカット2~5の<sup>2</sup>Dク ロマトグラムを示します。 カット2では、ピー クは検出されませんでした。 カット 3 は、 $^{1}D$  で はオーバーラップした化合物2と3が2つの 分離ピークとなることを示しています。カット 4は、大部分が化合物3ですが、ごく少量の化 合物2も含んでいます。カット5では、化合物 3のみが検出されました。



図 5.2 つの混合物、相対濃度 100 % の化合物 3 と相対濃度 0.05 % の化合物 1 と 2 (赤線)、 およびすべて相対濃度 0.05 % の 3 つの化合物 (青線) の <sup>1</sup>D クロマトグラム



図 6. 化合物 3 のピーク幅全体をサンプリングしたカット 1 ~ 9 の  $^{1}$ D クロマトグラム (上)。  $^{2}$ D クロマトグラム (下) はカット 2、3、4、5 の分析を示します。

ICH ガイドラインで定められた報告の必要な スレッシュホールドにおけるこのメソッドの有 用性を調べるために、3 つのクロロジフルオ ロ安息香酸異性体の各種混合物、相対濃度 100%の化合物3、相対濃度 0.05%、0.10%、 または 0.15%の化合物1 および2を調製しま した。図7に示すように、化合物3(3-クロロ -2,4-ジフルオロ安息香酸)および化合物2(5-クロロ-2,4-ジフルオロ安息香酸)のピーク面 積を Chemstationの2D-LC Viewer を用いて、す べてのカットの<sup>2</sup>D ピークの合計として計算し ました。表混合物は6回繰り返して分析しま した。表2に、3 つの異なる混合物について 計算したピーク面積の平均値と相対標準偏差 (RSD) を示します。低濃度の化合物2のピー クは、S/N比の平均値が混合物1、2、3でそれ ぞれ、3.2、5.9、12.2であり、これは定量下限 (LOO)以下でした。このことは、化合物2の積 分ピークが最大7.95%という比較的高いRSD 値からもわかります。化合物2は化合物3と 比較して相対面積比が0.03%、0.07%、0.14 %で検出されました。両方の化合物の<sup>1</sup>D分析 から求めたレスポンス係数が1.17であること を考慮すると、精度は71.6~108.5%と計算 できます。これは主要化合物と不純物の濃度 差が大きいことを考えると良好な結果といえ ます。このことは、一次元目でメインピークと 共溶出する不純物について、ICH ガイドライン 03A(R2) で報告が必要なスレッシュホールド (0.05 %) の相対濃度において高分離能サンプ リング 2D-LC を用いて API の不純物を分離お よび検出できることを示しています。



図 7. Chemstation の 2D-LC Viewer 画面。化合物 2 (5-クロロ-2,4-ジフルオロ安息香酸)、化合物 3 (3-クロロ-2,4-ジフルオロ安息香酸) のピーク面積を用いた クロロジフルオロ安息香酸の高分離能サンプリング 2D-LC 分析結果

表 2.3 つの異なる混合物を<sup>2</sup>D 検出器として光路長 10 mm の Max-Light カートリッジセルを備えた Agilent 1290 Infinity II DAD を用いて 6 回の連続分析の結果から算出したピーク面積の平均値と相対標準偏差。化合物 3 に対する化合物 2 の 面積比 (%)。<sup>1</sup>D ピークから算出したレスポンス係数を 1.17 として精度を計算しました。

3-クロロ-2,4- ジフルオロ安息香酸 (化合物 3)		5-クロロ-2,4- ジフルオロ安息香酸 (化合物 2)			面積		
	面積	RSD (%)	面積	RSD (%)	S/N	比 (%)	精度 (%)
混合物1(0.05:0.05:100%)	10,685.64	0.26	3.39	7.95	3.2	0.03	71.6
混合物 2 (0.10:0.10:100 %)	10,616.72	0.39	7.47	3.22	5.9	0.07	82.3
混合物 3 (0.15:0.15:100 %)	10,548.79	0.13	14.67	2.71	12.2	0.14	108.5

精度と真度をさらに向上させるために、 Agilent 1290 Infinity II HDR-DAD 不純物分析ソ リューションを <sup>2</sup>D 検出器として用いました。こ のソリューションでは、ダイナミック UV レンジ の直線範囲を広げるために、異なる光路長の Max-Light カートリッジセルを備えた 2 台のダ イオードアレイ検出器からのシグナルを組み 合わせます。高濃度の化合物 3 でも UV レン ジの直線範囲を超えることなく注入量を増や すことができました。S/N 比の平均値はすべて の混合物で LOQ を超えており、混合物 1、2、3 でそれぞれ 20.7、24.1、34.0 でした。化合物 2 の面積精度は大幅に向上し、RSD 値は 1.58 % 以下でした (表 3)。精度は 96.4 ~ 108.5 % で、 定量に最適です。

#### 脱アミドインスリンの分析

塩基性条件下でウシインスリンを脱アミド化 し、不純物を含む医薬品関連物質の実際のサ ンプルとしました。サンプルを HiRes サンプリン グ 2D-LC/MS で分析しました。Agilent OpenLab CDS ChemStation Edition と MassHunter からの



ることができます。  $\boxtimes$  8A に示すように、カット するシグナルを強調した <sup>1</sup>D UV クロマトグラム を MassHunter で表示することができます。  $\boxtimes$ 8B には MassHunter におけるフル 2D-MS を、  $\boxtimes$  8C にはフル 2D-UV シグナルを示します。10 のカットのそれぞれについて、MassHunter の <sup>2</sup>D Chromatogram Creator を用いて 1 つずつクロ

表3.3つの異なる混合物を<sup>2</sup>D 検出器として Agilent 1290 Infinity II HDR-DAD 不純物分析ソリューションを用いて 6回の連続分析の結果から算出したピーク面積の平均値と相対標準偏差。化合物3に対する化合物2の 面積比(%)。<sup>1</sup>D ピークから算出したレスポンス係数を1.17として精度を計算しました。

	3-クロロ-2,4- ジフルオロ安息香酸 (化合物 3)		5-クロロ-2,4- ジフルオロ安息香酸 (化合物 2)			面積	
	面積	RSD (%)	面積	RSD (%)	S/N	比 (%)	精度 (%)
混合物 1	50072.70	1.19	23.30	1.58	20.7	0.05	108.5
(0.05:0.05:100 %)							
混合物 2	49020.63	0.85	42.02	1.44	24.1	0.09	99.9
(0.10:0.10:100 %)							
混合物3	48698.74	1.34	60.37	1.47	34.0	0.12	96.4
(0.15:0.15:100 %)							



図 8. 脱アミドインスリン分析を Agilent MassHunter で表示したシグナル。A) HiRes サンプリングでの <sup>1</sup>D-UV シグナルのカット。B) フル <sup>2</sup>D-MS シグナル。C) フル <sup>2</sup>D-UV シグナル

マトグラムを作成しました。図9は、10のカットすべてについてのトータルイオンクロマトグ ラム (TIC) を重ねたものです。<sup>2</sup>D 分離の最初の 1 分間は、MS ダイバータバルブの切り替えに より廃液に流れるようにしました。この方法で フラクションに含まれる緩衝液塩が、MS イオ ン源に入らないようにしました。図9のそれぞ れの TIC において、分析時間1分でダイバー タバルブを切り替え、フローが MS に直接流れ るようになったことが視覚的にもわかります。 インスリンのメインピークが、カット2~10 に おいてリテンションタイム 2.06 分に異なる強 度で確認できます。 $^{2}D$  グラジエントには 90 %B でカラムを洗浄する短い時間があり、これによ り約 3.2 分にベースラインの急な増減が認め られます。図 10 にカット 7 の TIC と $^{2}D$ -UV シグ ナルの拡大図を示します。インスリンのメイン ピークは、一次元目からテーリングしていたこ とがわかります。 $^{2}D$  クロマトグラムでは、2 つ のピークが MS と UV シグナルの両方で確認で きます。図 11 に 2 つのピークの質量スペクト ルを抽出して示します。ピーク 1 の同位体比 は、インスリンの [M+5H]<sup>5+</sup> イオンとすることが できます。ピーク 2 の質量スペクトルは、イン スリンの単一の脱アミド生成物に相当する同 位体比 (中性分子質量のシフト 0.98 Da) を示し ています。同じ化合物がカット 6 でも検出され ました (データは非表示)。



図 9. Agilent MassHunter の <sup>2</sup>D Chromatogram Creator で作成したすべてのカットの <sup>2</sup>D クロマトグラム



図 10. カット 7 の <sup>2</sup>D クロマトグラム (MS (TIC) と UV シグナル (DAD2))



図 11. カット7の <sup>2</sup>D クロマトグラムにおけるピーク1と2の質量スペクトル

## 結論

このアプリケーションノートでは、高分離能サ ンプリング 2D-LC を用いて、構造的に類似して いる主要化合物のサンプルに含まれる低濃度 の不純物を検出できることを実証しました。ク ロロジフルオロ安息香酸の混合物を用いて、 新しい原薬の不純物についての ICH ガイドライ ンに従い、相対濃度が 0.05 %、0.10 %、0.15 %、 100%の<sup>1</sup>D分析で分離できない化合物を HiRes サンプリング 2D-LC により<sup>2</sup>D カラムに送 り、分離できることが示されています。5-クロ ロ-2,4-ジフルオロ安息香酸は 71.6~108.5% の精度で検出されました。面積精度は 2.71 ~ 7.95 % でした。<sup>2</sup>D 検出器として Agilent 1290 Infinity II HDR-DAD 不純物分析ソリューショ ンで注入量を増やした結果、精度が96.4~ 108.5%、面積精度が1.44~1.58%に向上し ました。脱アミドインスリンを実際のサンプル として、HiRes サンプリング 2D-LC/MS 分析をし ました。Agilent MassHunter の<sup>2</sup>D Chromatogram Creator を用いてデータを評価しました。インス リンのメインピークは一次元目ではテーリング していましたが、インスリンの単一の脱アミド 生成物が検出され、2つのカットにおける同位 体比で同定できました。

#### 参考文献

- Venkatramani, C. J.; *et al.* Assessing Stability-Indicating Methods for Coelution by Two-Dimensional Liquid Chromatography with Mass Spectrometric Detection. *J. Sep. Sci.* **2014**, *37*, 3214-3225.
- International Conference on Harmonization of Technical Requirements for Registration of Pharmaceuticals for Human Use – Impurities in New Drug Substances Q3A(R2), 2006.
- Petersson, P.; Haselmann, K.; Buckenmaier, S. Multiple Heart-Cutting Two Dimensional Liquid Chromatography Mass Spectrometry: Towards Real Time Determination of Related Impurities of Bio-Pharmaceuticals in Salt Based Separation Methods. J. Chromatogr. A 2016, 1468, 95-101.

- Stephan, S. High-Resolution Sampling 2D-LC with the Agilent 1290 Infinity II 2D-LC Solution. *Agilent Technologies Technical Overview*, publication number 5991-7637EN, 2016.
- Schneider, S. 2D-LC/MS Characterization of Charge Variants Using Ion Exchange and Reversed-Phase Chromatography Multiple Heart-Cutting 2D-LC Analysis of Innovator versus Biosimilar Monoclonal Antibodies. *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・テクノロジー株式会社 © Agilent Technologies, Inc. 2017 Printed in Japan, June 1, 2017 5991-8082JAJP

