バイオ医薬品

LC/UV および LC/MS による オリゴヌクレオチド分析のための 各種イオンペア試薬の評価

著者

Gerd Vanhoenacker, Cindy Lecluyse, Griet Debyser, Pat Sandra, Koen Sandra RIC Biologics Belgium

Sonja Schipperges, Sonja Schneider, Udo Huber Agilent Technologies, Inc. Germany

概要

オリゴヌクレオチド (ON) の LC/UV および LC/MS 分析に対する各種イオンペア試薬 (アミンおよび 酸性対イオン) の影響について検討しました。アミンの種類と対イオンの濃度が、ON のリテンションと 分解能に重要な役割を果たし、MS データに強く影響します。荷電状態と付加体の生成は、移動相の組 成の影響を受けます。本アプリケーションノートでは、ON の分析条件の最適化に多くの変数が利用で きることを示します。分析は、DAD および Agilent 6530 LC/Q-TOF による UV 検出を備えた Agilent 1290 Infinity II LC システムで実施しました。

はじめに

合成 ON は盛んに開発が進められており、さ まざまな疾患の治療に利用されています。 ON 治療薬は主に、核酸塩基、リボースまた はデオキシリボース糖(それぞれ RNA また は DNA)、およびリン酸からなる短い直鎖ポリ マーです。^{1,2} ON にはさまざまなサイズと形 のものがあり、治療効果を変更、強化し、特 定の標的に対する特異性を高め、安定性、デ リバリー、取り込みなどを改善するために、化 学修飾されることがよくあります。ON の種類 は幅広く、低分子干渉RNA (siRNA)、マイク ロ RNA (miRNA)、アンチセンスオリゴヌクレ オチド (ASO)、アプタマーなどのさまざまな クラスが生成されています。

ON は通常、ヌクレオチドを事前に定義された サイズ (つまり、ヌクレオチドの数 (n)) まで 1 つずつ追加することによって鎖を延長する多 段階プロセスで製造されます。合成はさまざ まな方法で行うことができ、一般に支持担体 と保護基の使用を伴います。ON 合成に関わ るすべてのステップで、不純物が発生する可能 性があります。よくある不純物には、反応が不 完全なショートマー (例えば、n-1、n-2) があ ります。サイクル数が増えるとターゲット ON の収量は低下します。あるいはロングマー(例 えば、n + 1) も形成され、脱アミノ化または 脱プリン化、酸化(典型的にはグアニン)、お よびその他の望ましくない反応から他の不純 物も牛じる可能性があります。製造時に混入 する不純物に加えて、ON の保管が分解生成 物につながる可能性があります。

主生成物の分析、純度の決定、不純物の定量 化・同定などの ON の分析は、化合物の性質 とサイズのために容易ではありません。近年、 合成 ON の種類と LC による分析に関する包 括的なレビューが発表されています。¹ ON は

本来、非常に極性の高い生成物であり、多く のアニオン基(通常、各ヌクレオチドに1つ のリン酸基を含む)を有しています。ON が修 飾されていない場合、または化学修飾による 極性全体への影響が小さい場合(一般的に あてはまります)、従来の逆相 LC では十分な リテンションが得られないため、他のモードを 適用する必要があります。アニオン交換クロマ トグラフィー (AEX)、親水性相互作用液体ク ロマトグラフィー (HILIC)、混合モードクロマ トグラフィー (MMC)、イオンペア液体クロマ トグラフィー (IPLC)、サイズ排除クロマトグラ フィー(SEC)など、さまざまな LC モードが ON の特性解析に適用されてきました。^{1,2} 上 記の組み合わせは、多次元で使用されること もあります。これは、包括的な 2D-LC³ または (複数の) ハートカット 2D-LC 設定で行うこ とができます。⁴ 要求される分析(QA/QC で あるか R & D であるか) に応じて、UV ベース の検出器または質量分析計を適用します。

ON 分析には、IPLC が群を抜いて最もよく使われる LC モードです。IPLC では、逆相固定相がイオンペア試薬を含む移動相と組み合わされます。これらの試薬は通常、アニオン性ON と相互作用して疎水性ペアを形成するアミンです。このペアはその後、逆相カラムに保持され、ON の長さ、種類、および化学修飾の有無に応じて分離されます。

移動相を適切に選択することが極めて重要 です。移動相は、アミン系イオンペア試薬と 酸性対イオンを決められた濃度と比率で正確 に、水中で混合することによって調製します。 この組み合わせにより、必要な ON の保持 が実現できます。メタノールやアセトニトリル などの有機溶媒を用いることで溶出が促進さ れ、グラジエント溶出により分離能の面で最 良の結果がもたらされます。特定のサンプル に対する結果は、選択した固定相、移動相添 加剤の濃度と種類、カラム温度、およびサン プル特性の複雑な相互作用によって決まりま す。クロマトグラフィーが移動相の組成によっ て影響を受けるだけでなく、LC/MS を使用す る場合、MS データ(感度、荷電状態、付加 イオンなど)も移動相の組成によって大きく影 響されます。

さまざまなグループが、ON のクロマトグラ フィーや質量分析の挙動に対する移動相の組 成の影響を調査してきました。⁵⁹ メソッドを注 意深く開発することで、分離能と感度の大幅 な改善が可能です。ただし、数多くの種類があ る ON に対して、最適に機能する一般的な単 一の条件セットを特定することは、不可能だと 考えられます。個々の ON に対して、イオンペ ア試薬の種類と濃度、対イオンの種類と濃度、 および双方の添加剤の濃度比を最適化するこ とにより、分離を最大化し、MS 検出を向上さ せることができます。

本アプリケーションノートでは、Agilent AdvanceBio オリゴヌクレオチドカラムを使用 し、さまざまな移動相で得られた結果を示し ます。分析は、ダイオードアレイ検出(DAD) および 6530 LC/Q-TOF を使用した 1290 Infinity II LC システムで実施しました。異な る移動相の組み合わせがクロマトグラフィー および質量分析の結果に与える影響を、DNA および RNA ON 参照サンプルおよび数種類 の治療用 ON により評価しました。データは、 ON 分析で従来より使用されてきた移動相に 代わる移動相があることを示しています。これ らの代替移動相は、同等以上の性能を示すだ けではなく、大幅に低いコストで分析を実施で きます。

実験方法

試料調製

次のアミンと酸は、Sigma-Aldrich(セン トルイス、ミズーリ州、米国)から購入しま した:トリエチルアミン(TEA)、n-ブチル アミン(BA)、ジブチルアミン(DBA)、n-ヘキシルアミン(HA)、酢酸トリエチルア ンモニウム(TEAA)、酢酸(AA)、および 1,1,1,3,3,3-ヘキサフルオロ-2-プロパノール (HFIP)。アセトニトリル(HPLC-S)、メタノー ル(ULC/MS)、および水(ULC/MS)は、 Biosolve(ファルケンスワールト、オランダ) から入手しました。サンプルは、ジエチルピ ロカーボネート(DEPC)処理水(Thermo Fisher Scientific、ウォルサム、マサチューセッ ツ州、米国)で希釈しました。

サンプル

- 標準混合物:
 - DNA ラダー標準 (部品番号 5190-9029)
 - RNA 分解能標準 (部品番号 5190-9028)
- サンプル:
 - 治療用 ON は、地元のバイオテクノロジー企業から入手しました。

移動相の前処理

酢酸移動相は、酢酸とそれぞれのアミンの等 モル溶液としました。500 mL を調製するた めに、酢酸を約 450 mL の水に加えて混合し ました。それぞれのアミンを加えて混合し、水 を加えて完全に 500 mL としました。この溶 液の pH は、少量の酢酸またはそれぞれのア ミンを添加して注意深<調整しました。

100 mM TEAA、 pH 7	酢酸:3 g トリエチルアミン:5.06 g		
100 mM HAA、 pH 7	酢酸:3 g ヘキシルアミン:5.06 g		
100 mM DBAA、 pH 7	酢酸:3 g ジブチルアミン:6.46 g		

HFIP 移動相は次のように調製しました。500 mL を調製するために、既定量の HFIP を約 400 mL の水に加えて混合しました。アミン をそれぞれ加えて混合し、水を加えて完全に 500 mL としました。これらの移動相はpH 調 整なしで使用しました。

	15 mM TEAA/400 mM HFIP pH 7.8 ~ 7.9	HFIP:33.6 g トリエチルアミン: 0.759 g
	15 mM TEAA/100 mM HFIP pH 8.6 ~ 8.7	HFIP:8.4 g トリエチルアミン: 0.759 g
	15 mM TEAA/25 mM HFIP pH 9.4 \sim 9.6	HFIP:2.1 g トリエチルアミン: 0.759 g
	15 mM HA/25 mM HFIP pH 9.4 ~ 9.6	HFIP:2.1 g ヘキシルアミン: 0.759 g
	15 mM DBA/25 mM HFIP pH 9.4 ~ 9.6	HFIP:2.1 g ジブチルアミン: 0.969 g

サンプル前処理

- DNA ラダー標準(部品番号 5190-9029)を DEPC 水1 mL に溶解しました。
- RNA 分解能標準(部品番号 5190-9028)を DEPC 水1 mL に溶解しました。
- サンプルを調製し、DEPC 水で希釈しました。

装置構成

Agilent 1290 Infinity II LC システムを Agilent 6530 LC/Q-TOF と組み合わせて使 用しました。

- 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)、サンプルサーモスタット (オプション 101)
- Agilent 1290 Infinity II マルチカラム サーモスタット(G7116B)
- Agilent 1290 Infinity II ダイオードアレイ 検出器(G7117B)、10 mm InfinityLab Max-Light カートリッジセル(G4212-60008)
- Agilent 6530 LC/Q-TOF (G6530A)、 Agilent Jet Stream ESI ソース

ソフトウェア

- LC-DAD 専用
 - Agilent OpenLab CDS ChemStation revision C.01.07 SR4 [505] 以降
- LC-DAD-MS 用
 - データ収集:機器コントロール用 Agilent MassHunter (B.08.00) 以降
 - データ解析: Agilent MassHunter Qualitative Analysis (B.07.00) 以降

一般的なメソッド設定

LC-DAD

パラメータ	設定値	
カラム	Agilent AdvanceBio オリゴヌクレオチド、2.1 × 50 mm、2.7 µm(部品番号 659750-702)	
移動相 A および B	結果と考察セクションを参照	
グラジエント	結果と考察セクションを参照	
流量	0.6 mL/min	
カラム温度	65 °C	
検出(DAD)	260/4 nm (リファレンス 355/20 nm) ピーク幅 > 0.025 min (10 Hz)	
注入 2 µL (ニードル洗浄フラッシュポート、3 秒間、メタノール)		
インジェクタ温度	12 °C	

LC(-DAD)/MS

パラメータ	設定値		
カラム	Agilent AdvanceBio オリゴヌクレオチド、2.1 × 50 mm、2.7 µm(部品番号 659750-702)		
移動相A	A 15 mM アミン/HFIP (25、100、または 400 mM) 水溶液		
移動相 B	メタノール		
グラジエント	結果と考察セクションを参照		
流量	0.6 mL/min		
カラム温度	65 °C		
検出 (DAD)	260/4 nm (リファレンス 355/20 nm) ピーク幅 > 0.025 min (10 Hz)		
注入	2 µL (ニードル洗浄フラッシュポート、3 秒間、メタノール)		
インジェクタ温度	6 °C		

LC/Q-TOF 設定

ソース設定	Agilent Jet Stream		
イオン化モード	ネガティブイオン化		
ドライガス温度	300 °C 8 L/min		
ドライガス流量			
ネブライザ圧力	35 psig		
シースガス温度	350 ℃		
シースガス流量	8 L/min		
ノズル電圧	1,000 V		
キャピラリー電圧	3,500 V		
取り込み設定			
フラグメンタ電圧	200 V		
質量範囲	400 ~ 3,200 m/z		
取り込みモデル	拡張ダイナミックレンジモード(2 GHz)		
スキャン速度	ン速度 3スペクトル/秒		
データ取り込み	セントロイドおよびプロファイル		

結果と考察

UV-DAD 検出用アミン-酢酸移動相

アミンー酢酸の組み合わせは、ON のLC/UV 分析に広く用いられている移動相添加物です。 一般に、100 mM 水溶液(アミンと酢酸の等 モル溶液)が使用され、分析成分をアセトニト リルによるなだらかなグラジエントを使用して 溶出させます。最も一般的な組み合わせは、 中性 pH の TEAA バッファ(トリエチルアミン (TEA) と酢酸)です。

TEAA は移動相添加物として一般的に受け入 れられていますが、特定の ON に対する性能 の点で必ずしも最良の選択肢であるとは限り ません。したがって、メソッド開発時には代替 のアミンを検討する必要があります。図 1 は、 さまざまな移動相を使用した DNA ラダーお よび RNA 分解能参照標準の LC/UV 分析を 比較したものです。TEAA 代替品のそれぞれ に対するグラジエントは、クロマトグラフィー 性能の直接比較を可能にするために、TEAA で得られたリテンションタイムと一致するよう にスケーリングしました。

酢酸と、ヘキシルアミン (HAA バッファ) ま たはジブチルアミン (DBAA バッファ) などの 代替アミンとの組み合わせは、参照サンプル の TEAA と比較して大幅に優れたパフォーマ ンスを示しています。図1の挿入図は分析の 短時間の詳細を示したものですが、分離効率 が高いことは明らかです。メインピークからの 不純物の検出と分離が大幅に改善されていま す。TEAA を使用する場合、DNA ラダー標準 は 10 ~ 14 % のアセトニトリルのグラジエン トでは 10 分間で溶出されますが、RNA 分離 標準はわずか6~8%のアセトニトリルで12 分間で溶出されます。RNA は、リボース糖に 2'-ヒドロキシルが存在するため、DNA に比べ てはるかに早く、またはかなり少ない有機溶 媒とともに溶出します。



図 1. IPLC 用のさまざまな移動相添加物を使用した 260 nm における LC/UV 分析。(A) DNA ラダー標準、(B) RNA 分解能標準。 移動相のグラジエントが図に示されています。その他のメソッドパラメーターについては、実験セクションを参照してください。

HAA および DBAA 移動相を使用した IPLC は、ON の保持がはるかに強いという特徴が あり、その結果、これらの条件下では有機溶 媒の含有量とグラジエントの勾配を増やさな ければなりません。性能向上はこれで本質的 に説明されます。これらのデータから、さまざ まな移動相添加物の評価をすることが、特定 の ON アプリケーションの分離を調整するの に有効なアプローチとなることは明らかです。

UV および MS 検出用アミン-HFIP 移動相

TEAA は IPLC/UV 分析による ON 分析には 有用ですが、ネガティブイオン化エレクトロ スプレーイオン化(ESI)MS 検出には適し ていません。20 年以上前に、Apffel ら^{5,6} が TEAA 移動相を使用した ESI-MS の感度不足 について報告しています。彼らは、HFIP がア セトニトリルに溶解しないため、有機溶媒とし てメタノールと組み合わせて HFIP を酸性対 イオンとして使用することを提案しました。こ の組み合わせが、ON の LC/MS 分析のリファ レンスとなっています。感度が高くなる主な根 本的な原因は、添加物の揮発性の違いにあ ります。HFIP は TEA よりも沸点が低く、エ レクトロスプレープロセス中に揮発性の高い HFIP が優先的に蒸発し、生成された液滴内 で枯渇します。その結果 pH が上昇してアルカ リ性となった条件下では、TEA-ON イオンペ アがソースで解離し、ON の検出能力が高まり ます。⁶ 酢酸は TEA よりも揮発性が低いため、 ESI 中の除去効率が低下します。その後、イオ ンペアの解離が少なくなるか、ソース内で安 定化することもあります。これは、MS の感度 を劇的に低下させます。

興味深いことに、UVと併用すると、TEAAと 比較してさらに優れたクロマトグラフィー性能 が TEA-HFIP の組み合わせで得ることができ ます。これは図 1 からも明らかです。図 1 で は、DNA ラダーと RNA 分解能標準の結果 を、以前に報告されたアミンとアセテートの組 み合わせと比較しています。15 mM TEA と 400 mM HFIP の組み合わせは、一般的に 組み合わせの開始点として使用されますが、 TEAA と比較して優れた性能を示し、DBAA と HAA をもとにした代替の移動相と同等の 性能を示します。TEA-HFIP の移動相のもう 1 つの利点は、TEAA と比較してシーケンス依 存性が少なく、ON のリテンションと溶出順序 が予測しやすいことです。⁷

ただし、HFIP を使用する本当の利点は、酢酸 を使用する場合と比較して、MS 感度が大幅 に向上することです。一方で HFIP の欠点は、 高品質の HFIP 試薬のコストが高いことです。 最適な MS 結果(すなわち、最高の感度と質 量スペクトル品質)を得るには、入手可能な 最高品質の HFIP を使用し、これらの移動相 を毎日調製することが推奨されます。HFIP を 大量に使用することを考えると(溶液を 400 mM 調製するのに約7g/100 mL が必要)、 これらの分析を定期的に実施するラボにとっ て、この消耗品のコストの増加はかなり大き なものになります。したがって、濃度が低い HFIP と移動相の組み合わせを検討すること は、分離の最適化のためだけでなく、分析あ たりのコストを削減するためにも、メソッド開 発の重要な部分となります。

UV および MS を用いて、アミンと HFIP のさ まざまな組み合わせを評価する実験を設定し ました。適用した移動相および、LC/UV スク リーニング実験中に使用したグラジエントの 概要を表1 にまとめます。前の実験とは対照 的に、メタノールのグラジエント勾配はすべて の分析で一定としました。15 mM TEA/400 mM HFIP の参照条件のリテンションウィンド ウに一致させる試みは行いませんでした。

表 1. アミンと HFIP の組み合わせとテストしたグラジエント

アミン(15 mM)	HFIP 濃度		
トリエチルアミン(TEA)			400 mM (=リファレンス)
ブチルアミン (BA)	25 mM	100 mM	
ジブチルアミン(DBA) [*]	2511111		評価せず
ヘキシルアミン(HA)			
pH バッファ	$9.45 \sim 9.55$	8.60 ~ 8.70	7.85
適用したグラジエント**			
グラジエント DNA ラダー標準お上びサンプル (3 % B/min)	10 ~ 55 % B	10~70 % B	10 ~ 55 % B
	15分	20 分	15分
グラジエント BNA 分解能標準 (2 % B/min)	$5\sim35$ % B	$5\sim45$ % B	$5\sim35$ % B
	15 分	20 分	15分

* アミン、特に DBA は、LC および MS ソースを汚染する傾向があります。 システムから汚染を除去するには、より広範なクリーン アップ手順が必要になる場合があります。

** 各グラジエントの最後に、洗浄ステップ(90% B、2分間)および再調整ステップ(最初の%B、5分間)を適用しました。

バッファ組成は、IPLC のリテンションに対して 大きな影響を及ぼします。図 2 に、テストを 行ったスクリーニングの組み合わせを用いて、 DNA ラダー標準の LC/UV 分析を実施した結 果を示します。ON のリテンションは HFIP 濃 度が低下すると一般に減少しますが、アルキ ルアミン鎖長が長くなるとリテンションは増加 します。RNA 分解能標準と治療用 ON につい ても同様の結果が得られました。アミン・酢酸 移動相スクリーニングから結論付けられるよ うに、アミン-HFIP を使用した移動相バッファ 組成の検討は、特定の ON アプリケーション の条件を最適化するために非常に効果的であ ることがわかります。 これらのスクリーニング結果に基づいて、移 動相の選択を LC/UV/MS を用いて評価しま した。デフォルト条件(15 mM TEA および 400 mM HFIP)および代替移動相で得られ た LC/MS クロマトグラムを図 3 ~ 6 に示し ます。グラジエントの開始条件と勾配は、特定 の移動相の組み合わせで分析したすべてのサ ンプルで一定としました。代替の移動相でリテ ンションが増加するので、次のステップではグ ラジエントを最適化して、特定のサンプルの分 離能、分析時間、および感度を向上させます。 これはここでは実施していません。 一般に、参照条件(15 mM TEA/400 mM HFIP)でのクロマトグラフィーは、代替の移動 相で行った結果と比較して優れたものではあ りません。リテンションは一般に、代替の移動 相(低濃度の HFIP)で増加します。これは、 分解能の向上に効果的です。参照条件と代替 条件に同じグラジエント勾配を使用した場合 に得られる分解能は、基本的に DNA ラダー 標準(図3)および RNA 分解能標準(図4) の影響を受けません。図3の質量スペクトル は、参照トリエチルアミン条件と比較して、代 替のへキシルアミン移動相では、DNA ラダー 標準の電荷エンベロープがより高い荷電種に シフトし、付加体が少ないことを示しています。







図 3. DNA ラダー標準の LC/MS 分析結果。15 mM TEA/400 mM HFIP 参照条件(左)および 15 mM HA/25 mM HFIP 代替移動相(右)。上: BPC、中:スペクトル 35mer、下:デコンボリュートしたスペクトル 35mer。移動相グラジエント:図に示されています。その他のメソッドパラメータ:実験セクションを参照してください。

RNA 分解能標準の分析では電荷状態のシフトは観察されませんでしたが、ここでも移動 相にジブチルアミンを使用する代替条件を使 用すると、付加体の少ないよりクリーンなスペ クトルが得られます。付加体の生成の減少は、 図3および図4の下部に示されているデコン ボリュートしたスペクトルで明瞭に示されてい ます。 治療用 ON サンプル A では、メインピークお よび、このメインピークの前に溶出する多くの 不純物ピークが示されています(図 5)。これ らのピークは、ショートマーに起因する可能性 があります。代替条件を使用すると、メインピー クと、メインピークのテールで溶出する不純物 (ロングマーではない)との間の分離度が向 上します。これは、HFIPの使用量が約8分の 1 で分解能が向上することを意味します。



図 4. RNA 分解能標準の LC/MS 分析結果。15 mM TEA/400 mM HFIP 参照条件(左)および 15 mM DBA/25 mM HFIP 代替移動相(右)。 上:BPC、中:スペクトル 21mer、下:デコンボリュートしたスペクトル 21mer。移動相グラジエント:図に示されています。 その他のメソッドパラメータ:実験セクションを参照してください。



図 5. サンプル A の LC/MS 分析結果。15 mM TEA/400 mM HFIP 参照条件(左) および 15 mM HA/25 mM HFIP 代替移動相(右)。 上:BPC、下:EIC メインおよびショートマー移動相グラジエント:図に示されています。その他のメソッドパラメータ:実験セクションを参照してください。 図 6 に示すのは、薬理学的に活性な 21mer および 22mer のアンチセンス鎖 (as) および 修飾された 21mer および 22mer のセンス 鎖(ss)で構成される 2 つの二本鎖 siRNA 治療薬(13.5~15.0 kDa、0.5 µM)の混合 物である、治療用 ON サンプル B の LC/MS データです。65 ℃の操作条件下では、二本鎖 が解離して一本鎖 (as)RNA と (ss)RNA が生 じることに注意してください。代替移動相(15 mM HA および 25 mM HFIP) を使用すると、 デフォルト条件(15 mM TEA および 400 mM HFIP) と比較して、リテンションが増加 し、21mer と 22mer (ss)RNA 間の分離能 が大幅に向上します。21 および 22mer (ss) RNA は、細胞への取り込みを実現するため に、リガンドで修飾されていることに注意して

ください。これらのクロマトグラフィーの観察 は、移動相とグラジエント条件が、その性質、 複雑さ、および分析の目的に応じて、任意の タイプの ON サンプルに対して最適化できるこ と、また最適化すべきであることを再度示して います。DNA ラダー標準について以前に観察 されたように、質量スペクトルは、代替条件お よびより少量の付加体で、電荷エンベロープ がより高い荷電種へシフトしていることを示し ています。

MS の結果は、HFIP の消費が少ないという 利点に加えて、一般的に代替移動相の方が 質の高い MS 結果をもたらすことを示してい ます。特に付加体の生成は、TEA-HFIP 参照 条件と比較して大幅に減少しています。付加 体の生成は、流路に存在する塩に対する ON

主鎖のリン酸ジエステルの親和性のためであ り、ON の IPLC/MS 分析でよく知られた現象 です。付加体(主にナトリウムとカリウム)の 量は、いくつかの要因の影響を受けます。移 動相、特に HFIP の品質は、付加体の有無に 大きな影響を与えます。このため、移動相は 新たに調製して使用し、入手可能な最高品質 の HFIP を使用することが推奨されます。本 実験では 純度 99 % 以上の HFIP を使用しま した。より高い純度のものが入手可能ですが、 それに応じてコストが増加します(約5倍)。 移動相のボトルや LC システムとキャピラリは すべて、できる限り清浄にする必要がありま す。MS スペクトルの品質を向上させ、付加体 の量を減らすために、LC システムの前処理ま たは洗浄が必要になる場合があります。



図 6. サンプル B の LC/MS 分析結果。15 mM TEA/400 mM HFIP 参照条件(左)および 15 mM HA/25 mM HFIP 代替移動相(右)。 上:BPC、中:スペクトル (as) RNA 21mer および 22mer、下:デコンボリュートしたスペクトル (as) RNA 21mer および 22mer。 移動相グラジエント:図に示されています。その他のメソッドパラメータ:実験セクションを参照してください。

結論

さまざまな ON の分析において、移動相の選 択がクロマトグラフィーと質量分析の両方の 挙動に重要な役割を果たすことを示す結果が 得られました。従来の移動相の組成は、一般 的な ON 分析でその価値が実証されています が、移動相の組成は特定のサンプルに対して 最適化できることが実証されました。LC/UV に一般的に使用される酢酸ベースの移動相 は、主にトリエチルアミンと酢酸の組み合わせ に基づいています。本アプリケーションノート のデータは、クロマトグラフィーの品質を向上 させるには多くの場合、ヘキシルアミンやジブ チルアミンなどの代替アミンを使用する方が 優れていることを示しました。LC/MS の作業 では、感度が向上するためHFIP ベースの移 動相が推奨されます。ここでも、アミンの種類 の選択と HFIP の濃度が結果に大きく影響し ます。移動相中の高価な HFIP の濃度を減ら すことのもう 1 つの利点は、日常的に ON を 分析する場合の分析コストの削減です。

参考文献

- Goyon, A.; Yehl, P.; Zhang, K. Characterization of Therapeutic Oligonucleotides by Liquid Chromatography. J. Pharma. Biomed. Anal. 2020, 182, 113105.
- Vanhinsbergh, C. J. Analytical Separation Methods for Therapeutic Oligonucleotides. LC-GC, Chromatography Online, Advances in Biopharmaceutical Analysis (Special Issue) 2020, 33(10), 20–26.
- Li, et al. Comprehensive Hydrophilic Interaction and Ion-Pair Reversed-Phase Liquid Chromatography for Analysis of Di- to Deca-Oligonucleotides. J. Chromatogr. A 2012, 1255, 237-243.
- Krieger, S.; Dickhut, C. Direct Analysis of In-Process Oligonucleotides Without Manual Purification. *Agilent Technologies application note*, publication number 5991-9490EN, **2018**.
- Apffel, A. et al. Analysis of Oligonucleotides by HPLC-Electrospray Ionization Mass Spectrometry. Anal. Chem. 1997, 69, 1320-1325.
- Apffel, A. et al.New Procedure for the Use of High-Performance Liquid Chromatography-Electrospray Ionization Mass Spectrometry for the Analysis of Nucleotides and Oligonucleotides.J. Chromatogr.A 1997, 777, 3-21.

- Gilar, M. *et al*.lon-Pair Reversed-Phase High-Performance Liquid Chromatography Analysis of Oligonucleotides: Retention Prediction.*J. Chromatogr.A* 2002, 958(1-2), 167-182.
- McGinnis, A. C.; Grubb, E. C.; Bartlett, M. G. Systematic Optimization of Ion - Pairing Agents and Hexafluoroisopropanol for Enhanced Electrospray Ionization Mass Spectrometry of Oligonucleotides.*Rapid Commun.Mass* Spectrom.**2013**, 27(23), 2655–2664.
- Gong, L.; McCullagh, J. S. O. Comparing Ion-Pairing Reagents and Sample Dissolution Solvents for Ion-Pairing Reversed-Phase Liquid Chromatography/Electrospray Ionization Mass Spectrometry Analysis of Oligonucleotides.*Rapid Commun.Mass Spectrom.***2014**, 28(4), 339-350.

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. 2021 Printed in Japan, January 4, 2021 5994-2957JAJP DE44180.3438773148

