

高分解能 LC/MS を用いた mRNA ポリ A 配列変異体の分析

著者

Brian Liau Agilent Technologies, Inc.

はじめに

2020 年の SARS-CoV-2 パンデミックにより生じた緊急事態の際には、政策立案者も製薬会社も同様 に mRNA ワクチンを前例のない速さで開発して配布することが要求されました。mRNA ワクチンは臨 床試験では優れた安全性と有効性を示しており^{1~4}、代替技術を基礎にしたワクチンよりも性能が優れ ています。mRNA ワクチンは遺伝子治療効果があると考えられているため⁵、FDAのガイダンスでは製 品関連の不純物を幅広く特性解析することが求められています。この中には、配列にわずかな異常があ る mRNA 分子の個体群が含まれている場合があり、これは配列変異体として知られています。さらに、 mRNA ワクチンでは最適な安定性と生物活性を実現するために、3' 末端において A ヌクレオチドの非 常に長い反復部分 (ポリ A) が必要になります。⁶そのため、ポリ A 配列の長さと組成はともに重要品 質特性です。

本研究では Agilent AdvanceBio 6545XT LC/Q-TOF を使用して、in vitro 転写系の一 般的な成分である大腸菌ポリ A ポリメラーゼ (PAP) によって形成されるポリA テール配 列を分析します。分析結果から、PAP は ATP に対して完全には選択的ではなく、CTP およ び UTP プリカーサの両方に影響を及ぼすこ とにより、標準の in vitro 転写条件下で大量の 望ましくない C および U ヌクレオチドを取り 込むことがわかっています。これらの配列変異 体は製品関連の不純物と見なすことができる ため、分析結果は PAP の使用に対して警告 を与えています。また LC/MS は核酸治療の プロセス最適化と品質管理において、高感度 で効率的な評価の高いメソッドであることを示 しています。

使用した略語:

- ATP アデノシン三リン酸
- CTP シチジン三リン酸
- UTP ウリジン三リン酸
- GTP グアノシン三リン酸
- A、C、U、および G ヌクレオチド アデノ
 シン、シチジン、ウリジン、およびグアノシン
 ンーリン酸
- ポリA-ポリアデノシン
- PAP 大腸菌ポリ A ポリメラーゼ
- RNA-seq-RNA シーケンシング

実験方法

mRNA の in-vitro 転写

上流 T7 プロモーターおよび下流 BGH ター ミネーター配列が側面に配置されている 3822 nt 遺伝子をコード化する pCMV3 プ ラスミドは Sino Biological から購入しまし た。DNA 配列は、T7 および BGH ターミネー タープライマー (Agilent ヘラクレス II Fusion DNA ポリメラーゼ、p/n 600677)を使用し て、35 サイクルで PCR 増幅しました。 クリー ンアップ後 (Agilent StrataPrep PCR 精製 キット、p/n 400771)、増幅した dsDNA を Agilent 2100 バイオアナライザと Agilent DNA 7500 キット (p/n 5067-1506) で分 析して濃度を測定し、増幅の均一性を評価し ました。次に、 増幅した dsDNA (~13 nM) を HiScribe T7 ARCA mRNA キット (New England Biolabs、p/n E2060S)を使用し て in vitro 転写し、メーカー推奨のプロトコル によって含有 PAP 酵素でテーリングしてから、 LiCl 沈殿させました。PAP テーリングの前後 に、転写 mRNA を Agilent 2100 バイオアナ ライザと Agilent RNA 6000 ナノキット (p/ n 5064-1511) で分析し、反応をモニタリング しました。

図 1A に示されているように、PAP の選 択性実験では 5' および 3'-OH を含む合 成 10-mer ポリA 配列 (Integrated DNA Technologies) を、反応ごとにただ 1 つの プリカーサヌクレオシド三リン酸 (1 mM の ATP、CTP、UTP、GTP のいずれか)を使用 して、37 °Cで 30 分間 PAP 酵素により拡張し ました。

サンプル前処理

約 20 ピコモルのポリ A テーリング mRNA を 37 °Cで 3 時間 1,000 U の RNase T1 で 消化し、ポリ A 配列を遊離しました。各サン プルを 200 µL のオリゴ-dT 磁気ビーズを使 用して 5 回精製し、ポリ A 配列をプルダウン しました。⁷各プルダウンを 50 µL の 1x IDTE バッファ(Integrated DNA Technologies、 p/n 11-05-01-05)に 溶 出 し、 最 終 量 250 µL にプールしました。LC/MS 分析の 前に、プールした溶出液を Vivaspin 500 カートリッジと 10 kDa MWCO(Sartorius、 p/n VS0102)を使用して 60 µL の脱イオン 水に脱塩しました。



Poly-A tailing experiments



図 1. このアプリケーションノートで実施したテーリング反応の概略図。(A)反応ごとに1つのプリカーサのみを 使用して実施した RNA プライマーでの反応。(B)標準条件下において *in vitro* 転写 mRNA で実施した反応 (すべてのプリカーサ)

ポリ A 配列の LC-DAD/MS

機器の構成は次のとおりです。

- 1290 Infinity II LC とダイオードアレイ検 出器 (p/n G7117B)
- 6545XT AdvanceBio LC/Q-TOF

流路から注意してガラスを取り除き、アルカリ 金属付加物を減少させました。移動相の容器 として Agilent ナルゲンボトル (p/n 9301-6460)を使用し、各溶媒ラインにステンレスフ リットを備え付けました。Agilent ポリプロピレ ンサンプルバイアル (p/n 5190-2242)を使 用しました。最初に使用する前に、LC システ ムとカラムに 50 % MeOH + 0.1 % ギ酸溶液 を一晩通液し、アルカリ金属付加物をさらに減 少させました。必要な場合に、実験間でシステ ムを清掃するには通常、50 % MeOH + 0.1 % ギ酸による 30 分間の通液で十分でした。⁸

ポリA 配列を PLRP-S カラム (2.1 × 50 mm、5 μ m、1,000 Å、p/n PL1912-1502) で分離しました。クロマトグラフィー分離能を向上させるために、PAP 選択性実験ではAgilent Infinity Poroshell 120 HPH-C18 カラム (2.1 × 50 mm、1.9 μ m、120 Å、p/n 699675-702)を使用しました。移動相と LC グラジエントを表 1 に示します。質量分析計をネガティブイオンモードで表 2 の設定にして操作し、BioConfirm 10.0 を表 3 のデコンボリューション設定にしてデータを解析しました。

表1.移動相とLC グラジエントの設定

Agilent 1290 Infinity II LC システム							
カラム	Agilent InfinityLab Poroshell 120 HPH-C18、 1.9 µm、2.1 × 50 mm、120 Å	Agilent PLRP-S、5 μm、2.1 × 50 mm、 1,000 Å					
溶媒 A	15 mM ジブチルアミン + 25 mM HFIP 脱イオン水溶液						
溶媒 B	15 mM ジブチルアミン + 25 mM HFIP メタノール溶液						
グラジエント	0~2分で15%B 12分で30%B 12.1~13分で90%B	0~1分で15%B 10.5分で45%B 10.6~11.5分で90%B					
カラム温度	50 °C	80 °C					
流量	0.4 mL/min						
注入量	10 ~ 20 μL						

表2. 質量分析計の設定

Agilent 6545XT AdvanceBio LC/Q-TOF								
	LC/MS	LC/MS/MS						
取り込みモード	ネガティブ、標準(3,200 m/z)質量範囲、高感度(2 GHz)							
ガス温度	350 °C							
ガス流量		12 L/min						
ネブライザ		55 psig						
シースガス温度		275 °C						
シースガス流量		10 L/min						
Vcap		4,500 V						
ノズル電圧		2,000 V						
フラグメンタ	250 V							
スキマ電圧	65 V							
MS1 範囲	400 ~ 3,200 m/z							
MS1 スキャン速度	2 Hz	5 Hz						
MS2 範囲		50 ~ 3,200 m/z						
MS2 スキャン速度		3 Hz						
MS2 選択幅:		中 (~ 4 amu)						
コリジョンエネルギー		0、40、60 V						
MS2 のスレッシュホールド		オン、3 回繰り返して 0.2 分間排除						
プリカーサアバンダンスベースの	_	あり						
スキャンスピード								
ターゲット(カウント/スペクトル)		25,000						
MS2 累積時間制限の使用		あり						
純度		100 % ストリンジェンシー、30 % カットオフ						
プリカーサの順位		+3、+2、+1 のみのアバンダンスによる						
リファレンス質量		1,033.9881						

表3. デコンボリューションの設定

Agilent MassHunter BioConfirm B10.0 の設定						
オリゴヌクレオチドの長さ	≤ 30 nt	≥ 90 nt				
デコンボリューションアルゴリズム	最大エントロピー					
ベースライン減算	1					
付加物	プロトン損失					
質量範囲	3,000 ~ 10,000 Da	$30,000 \sim 60,000$ Da				
質量ステップ	0.05 Da	0.05 Da				
限られた m/z 範囲の使用	1,040 ~ 3,200	800~2,500				

結果と考察

最初の試験では、1 mM の ATP が存在する 中で 10 個の反復 A ヌクレオチド (A_{10}) で構 成される合成 RNA プライマーで、PAP によ り拡張されたポリ A 配列を分析しました。図 2 に示すように、これにより二峰性分布のポリ A 配列が生成されましたが、1 つの個体群は 2.5 ~ 6 分に溶出する短鎖オリゴヌクレオチ ドで構成されており、もう1つの個体群は~ 10.6 分に広いピークで溶出する長鎖ヌクレオ チドで構成されていました。質量分析では、サ イズが11~22 ntの範囲の短鎖個体群を示 した(図3D)一方で、長鎖個体群は長さが 108~149 ntの範囲でした(35,492.89~ 48,990.17 Da、図4C)。

短鎖オリゴヌクレオチド個体群で選択した3 つのピークから抽出してデコンボリュートした 質量スペクトルを図3に示します。質量スペク トルは、主にプロトン損失から生成された二 価イオンと三価イオン、およびナトリウム付加 物の少数個体群で構成されていました。同位 体分離してデコンボリュートした質量スペクト ルは、それぞれのモノアイソトピックピークに 基づいて誤差 < 5 ppm により A₂₀、A₂₁、およ び A₂₂ と同定されました(図3B~3D)。



図 2. ATP のみが存在する中での 260 nm の UV 吸光度(A:リファレンス = 360 nm)および PAP で拡張された RNA プライマーのトータルイオンクロマトグラム(B)。 分離は PLRP-S カラムで実施しました。



図 3. PAP によって形成された A₁₁ ~ A₂₂ オリゴヌクレオチド。(A) 選択した 3 つのピークと抽出された質量スペクトルを示す (B ~ D) トータルイオンカレントクロマトグラム。 A₂₀ (M_{obs} = 6,519.12 Da、M_{theo} = 6,519.09 Da)、A₂₁ (M_{obs} = 6,848.16 Da、M_{theo} = 6,848.15 Da)、A₂₂ (M_{obs} = 7,177.22 Da、M_{theo} = 7,177.20 Da) のデコンボリュート した質量スペクトルを挿入図として示しています。M_{obs}: 観察値のモノアイソトピック質量、M_{theo}: 理論上のモノアイソトピック質量

長鎖オリゴヌクレオチドの一部をサンプリング してデコンボリューションしました(図 4A)。 10~10.3分に抽出された質量スペクトル の電荷エンベロープは主に800~2,500 m/zに収まっており(図 4B)、目的の質量範 囲 30~60kDaにデコンボリュートしまし た。デコンボリュートした質量スペクトル(図 4C)は明らかに、34~50kDaのサンプル ピークの不均一な個体群を示していました が、329.2±1Daごとに均一に分離されまし た(図 4D)。この質量増分はAヌクレオチド の単一添加の場合と一致しており、329.209 Daごとに理論上の平均質量が増加していま す。表 4 では、図 4D の質量ピークは、理論 上の質量と観察値の質量が ≤ 1.16 Da の差で A₁₂₁ ~ A₁₃₈ として高い信頼性でアノテーショ ンされたことを示しています。

ATP に対する PAP の選択性を評価するため に、1mM の CTP、UTP、または GTP のみが 存在する中で PAP を RNA プライマーに加え て同じ実験を実施しました。長ポリマー鎖の 拡張は観察されませんでしたが、C ヌクレオ チドの最大 2 個の単量体 (図 5A) または 1 個の U ヌクレオチド (図 5B) の RNA プライ マーへの添加をクロマトグラフ的に分離してお り、これは ATP に対して PAP が完全には選 択的ではなかったことを示しています。この実 験ではグアノシンーリン酸の添加は観察され ませんでしたが(図 5C)、長い反応時間また は高 GTP 濃度において感知できる量が添加 された可能性は除外できませんでした。全体 的に、PAP は ATP で最高の活性を示してお り、続いて降順に CTP、UTP、GTP でした。



図 4. PAP によって形成された A₁₀₈ ~ A₁₄₉ オリゴヌクレオチド。(A) デコンボリューションのためにサンプリングされた領域 10 ~ 10.3 分を示すトータルイオンカレント クロマトグラム。サンプリング領域の(B)電荷エンベロープと(C) デコンボリュートした質量スペクトル。破線の矢印(左側 = 35,492.89 Da、右側 = 48,990.17 Da)は、 高い信頼性で A₁₀₈ ~ A₁₄₉ と同定できた質量ピークの範囲を示しています。(D) 39,773.03 ~ 44,710.28 Da のピーク間で一定の間隔を示す拡大してデコンボリュートした 質量スペクトル

表	4.	义	4D	から	5ア.	ノテ・	ーショ	ョン	され	った	質量	ピー	ク
---	----	---	----	----	-----	-----	-----	----	----	----	----	----	---

オリゴヌクレオチド	観察値の質量(Da)	理論上の質量 (Da)	質量差(Da)
A ₁₂₁	39,773.03	39,772.08	0.95
A ₁₂₂	40,101.53	40,101.28	0.25
A ₁₂₃	40,431.09	40,430.49	0.6
A ₁₂₄	40,759.75	40,759.70	0.05
A ₁₂₅	41,089.46	41,088.90	0.56
A ₁₂₆	41,418.82	41,418.11	0.71
A ₁₂₇	41,748.39	41,747.32	1.07
A ₁₂₈	42,077.06	42,076.52	0.54
A ₁₂₉	42,406.35	42,405.73	0.62
A ₁₃₀	42,735.87	42,734.94	0.93
A ₁₃₁	43,065.04	43,064.15	0.89
A ₁₃₂	43,393.78	43,393.35	0.43
A ₁₃₃	43,723.09	43,722.56	0.53
A ₁₃₄	44,052.93	44,051.77	1.16
A ₁₃₅	44,381.54	44,380.97	0.57
A ₁₃₆	44,710.28	44,710.18	0.1



図 5. (A) GTP、(B) UTP、(C) CTP に対する PAP の無差別な活性を示す UV 吸光度 (Abs = 260 nm、Ref = 360 nm) クロマトグラム。グアノシンーリン酸の添加物 は検出されませんでした。 (D) パネル (C) に示されている A₁₀、A₁₀C、および A₁₀CC の相対的な定量。分離は Agilent Poroshell 120 HPH-C18 カラムで実施しました。

未修飾 RNA プライマーおよび C または U ヌ クレオチドで拡張してデコンボリュートした質 量スペクトルを図 6 に示します。同位体分離 してデコンボリュートした質量スペクトルは、 それぞれのモノアイソトピックピークに基づい て誤差 < 13 ppm により A_{10} C、 A_{10} CC、 および A_{10} U と同定されました。MS/MS 実 験では、C および U ヌクレオチドを実際に RNA プライマーの 3' 末端に加えると(図 7)、特徴的な二価 y-イオン 1601.271 m/z と 1601.758 m/z を形成したことが示されまし た。一方、未修飾 RNA プライマーを 3' A ヌク レオチドで末端化すると、フラグメンテーショ ンで二価 y-イオン 1448.749 m/z を生成しま した。 次に、RNA 6000 ナノキットを搭載したバイオ アナライザと LC/MS により、全長、in vitro 転 写 mRNA を分析しました。テーリングの前、 転写 mRNA は予測した長さ~3,800 nt を示 し、PAP との反応後~4,200 nt まで増大し ましたが(図 8A)、これはポリ A のテーリン グに成功したことを示しています。



図 6. (A) 未修飾 A₁₀ RNA プライマー (M_{obs} = 3,228.61 Da、M_{theo} = 3,228.57 Da)、(B) 1 個の C ヌクレオチドで拡張 (M_{obs} = 3,533.65 Da、M_{theo} = 3,533.61 Da)、 (C) 2 個の C ヌクレオチドで拡張 (M_{obs} = 3,838.67 Da、M_{theo} = 3,838.65 Da)、(D) 1 個の U ヌクレオチドで拡張 (M_{obs} = 3,534.62 Da、M_{theo} = 3,534.59 Da)の 抽出してデコンボリュートした質量スペクトル。M_{obs}: 観察値のモノアイソトピック質量、M_{theo}: 理論上のモノアイソトピック質量



図7. 異なる3'末端の特徴である診断イオンを示す選択したオリゴヌクレオチドの MS/MS

全長 mRNA サンプルを RNase T1 で消化し、 次にオリゴ dT 磁気ビーズで繰り返しプルダウ ンして精製テール配列を生成しました。PAP 拡張 RNA プライマーと同様に、*in vitro* 転写 mRNA から生成されたテール配列は、3.7 ~ 7.5 分に溶出するオリゴヌクレオチドの短鎖 個体群と~10.6 分に溶出する長鎖個体群の 両方で構成されていました(図 8B)。短鎖個 体群で選択したピークから抽出してデコンボ リュートした質量スペクトルは、長さが 16 ~ 27 nt の範囲のポリ A 配列を示しており、そ れぞれに誤って取り込まれた単一の U ヌクレ オチドが含まれていました (図 9)。このデー タセットでは観察されていませんが、他の実験 でも誤って取り込まれた C ヌクレオチドが観 察されました。

M. Beverly et al.⁷ で示されているように、PAP によって形成されたテール配列は、遺伝学的 に鋳型化されているポリA 配列と比較して長 さがかなり不均一です。分析結果から、この 不均一性は 4 種類のプリカーサヌクレオシド 三リン酸がすべて存在する標準条件下でテー リング反応が発生した際に、異なる数の C お よび U ヌクレオチドが誤って取り込まれたこと によって生じており、長鎖テール配列の質量ス ペクトルをデコンボリュートするのは非常に困 難であることがわかります。



図8. (A) PAP でテーリング前(レーン1)およびテーリング後(レーン2)の *in vitro* 転写 mRNA のバイオアナライザ分析。(B) 4 種類のヌクレオシドリン酸プリカーサが すべて存在する中での 260 nm の UV 吸光度(上部パネル)および *in vitro* 転写 mRNA に追加されたポリ A 配列のトータルイオンクロマトグラム(下部パネル)。 分離は PLRP-S カラムで実施しました。



図 9. U ヌクレオチドが誤って取り込まれた A₁₆ ~ A₂₇ オリゴヌクレオチド。(A) 選択した 3 つのピークと抽出された質量スペクトルを示す (B ~ D) トータルイオンカレント クロマトグラム。A₁₈ + U (M_{obs} = 6,167.04 Da、M_{theo} = 6,167.01 Da)、A₁₉ + U (M_{obs} = 6,496.10 Da、M_{theo} = 6,496.07 Da)、A₂₀ + U (M_{obs} = 6,825.15 Da、 M_{theo} = 6,825.12 Da) のデコンボリュートした質量スペクトルを挿入図として示しています。M_{obs}: 観察値のモノアイソトピック質量、M_{theo}: 理論上のモノアイソトピック質量

結論

本研究で次のことが示されました。(1) 長鎖 (121 ~ 136 nt) 不均一ポリA 配列のイン タクト質量は、質量スペクトル全体をデコンボ リューションすることによって正確に測定でき ます。(2) PAP は標準の *in vitro* 転写条件下 で ATP に対して完全には選択的ではなく、C および U ヌクレオチドの両方がポリA テール 配列に加えられています。 in vitro 研究ではこれらの配列変異体は重要 でない場合がありますが、規制当局の観点か らすると非常に重要です。特に、T7 ポリメラー ゼのような他の in vitro 転写酵素は、ずれや転 写停止のようなメカニズムを通して配列変異 体を生成する場合もあり⁹、これらの不純物を 検出するための高感度で選択的なメソッドが 必要であることを明確に示しています。 このような高感度と選択性を実現するために、 過去のある研究では放射能標識ヌクレオチド を使用した PAP のオフターゲット活性につい て示しました。¹⁰製造環境では、このような手 法は危険で不適切な場合があります。LC/MS は、このような試薬を使用しなくても単一ヌ クレオチドの選択性を実現できます。さらに LC/MS は、バイアスやアーティファクトをもた らすことで知られる、RNA-seq の特徴である 非常に長い逆転写、ライゲーション、および増 幅のステップなしに、配列変異体を検出して 定量できます。

参考文献

- Jackson, L. A. et al. An MRNA Vaccine against SARS-CoV-2 — Preliminary Report. N. Engl. J. Med. 2020, 383, 1920–1931.
- Mulligan, M. J. et al. Phase I/II Study of COVID-19 RNA Vaccine BNT162b1 in Adults. *Nature* 2020, 586, 589–593.
- Pfizer and BioNTech Announce Vaccine Candidate Against COVID-19 Achieved Success in First Interim Analysis from Phase 3 Study.
- 4. Promising Interim Results from Clinical Trial of NIH-Moderna COVID-19 Vaccine.
- Wadhwa, A. *et al.* Opportunities and Challenges in the Delivery of MRNA-Based Vaccines. *Pharmaceutics* **2020**, *12*, 102.
- Kahvejian, A. *et al.* Mammalian Poly (A)-Binding Protein Is a Eukaryotic Translation Initiation Factor, Which Acts via Multiple Mechanisms. *Genes Dev.*2005, *19*, 104–113.

- Beverly, M.; Hagen, C.; Slack, O. Poly A Tail Length Analysis of *In Vitro* Transcribed MRNA by LC/MS. *Anal. Bioanal.Chem.* 2018, 410, 1667–1677.
- Birdsall, R. E. *et al.* Reduction of Metal Adducts in Oligonucleotide Mass Spectra in Ion-Pair Reversed-Phase Chromatography/Mass Spectrometry Analysis. *Rapid Commun. Mass Spectrom.***2016**, 30, 1667–1679.
- 9. Tateishi-Karimata, H.; Isono, N.; Sugimoto, N. New Insights into Transcription Fidelity: Thermal Stability of Non-Canonical Structures in Template DNA Regulates Transcriptional Arrest, Pause, and Slippage. *PLOS ONE* **2014**, *9*.
- Yehudai-Resheff, S.; Schuster, G. Characterization of the E.Coli Poly(A) Polymerase: Nucleotide Specificity, RNA-Binding Affinities and RNA Structure Dependence. *Nucleic Acids Res.* 2000, *28*, 1139–1144.

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, March 8, 2021 5994-3005JAJP RA44231.5456365741

