

高分解能 LC/MS を用いた mRNA 5['] キャッピングの高速分析

はじめに

SARS-CoV-2 の大流行により、mRNA の工業的な生産に注目が集まっています。mRNA ワクチンはウ イルスに対して最も効果的な手段の 1 つであることが証明されており、*in vitro* 転写により合成された 高品質の mRNA に依存しています。¹mRNA がタンパク質に効率的に翻訳されるかどうかは、転写プ ロセス中²または転写プロセス後³に mRNA に付加されるキャッピングと呼ばれる 5'-末端ジヌクレオチ ド修飾に大きく依存しています。ここで重要なのは、正常にキャッピングされる物質の割合と付加された キャッピング構造の種類は、インプットする物質の品質、反応条件、mRNA 配列などの要因に影響され るということです。このため mRNA 5' キャッピングは重要な品質特性であり、十分な特性解析とモニタ リングが必要です。

今回の実験では Agilent 6545XT AdvanceBio LC/Q-TOF を使用して、mRNA キャッピングの高速分 析とサンプル処理を組み合わせたところ、分析時間はわずか 75 分でした。高反応温度で耐熱性酵素を 使用して、キャッピング構造を含む 5' 末端オリゴヌクレオチドを遊離することにより、サンプル前処理を 高速化しています (図 1)。次に、シンプルなクリーンアップ手順のみを使用して、キャッピングされたオ リゴヌクレオチドを AdvanceBio オリゴヌクレオチドカラムでサンプルマトリックスから分離します。

このメソッドの有用性を実証するために、分析困難な mRNA 配列に対してワクシニア酵素キャッピング 反応を最適化しました。この結果は、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 的なメソッドとして高分解能 LC/MS が有用であることを示しています。

使用されている略語

- DNA:デオキシリボース核酸(Deoxyribose nucleic acid)
- PCR:ポリメラーゼ連鎖反応 (Polymerase chain reaction)
- mRNA:メッセンジャーリボ核酸(Messenger ribonucleic acid)
- ARCA:抗リバースキャップアナログ (Anti-reverse cap analogue)
- UTR:非翻訳領域(Untranslated region)
- GTP: グアノシン三リン酸(Guanosine triphosphate)
- NTP: ヌクレオシド三リン酸 (Nucleoside triphosphate)
- SAM: S-アデノシルメチオニン (S-adenosylmethionine)

著者

Brian Liau Agilent Technologies, Inc.

実験方法

PCR プライマーと合成 RNA/DNA キメラプローブはすべて、Integrated DNA Technologies で合成しました。

In vitro 転写と mRNA キャッピング

T7 プロモーター下流のモデル~3,800 nt 配列をコード化するプラスミド は Sino Biological から購入しました。配列には、5' および 3' UTR とコー ディング領域が含まれていました。目的の T7 プロモーターと配列は、 ヘラクレス II Fusion DNA ポリメラーゼキット(Agilent p/n 600677) を使用して PCR 増幅し、StrataPrep PCR 精製キット(Agilent p/n 400771)を使用してクリーンアップしました。PCR 反応の特異性と 生成物の濃度を、Agilent 2100 バイオアナライザ機器と Agilent DNA 7500 キット(p/n 5067-1506)で測定しました。

次に、メーカー推奨のプロトコルで T7 RNA ポリメラーゼ (New England Biolabs M0251) を使用して、PCR 生成物を mRNA に転写しました。100 µL *in vitro* 転写反応物を 100 µL の nuclease-free water で希釈した後、DNase-I (New England Biolabs M0303) 分解により DNA テンプレートと残留 PCR プライマーを除去しました。次に、70 µL の 8 M LiCl を加えて mRNA を沈殿させ、 $-20 \degree C$ で一晩冷却してから、12,000x g で 15 分間 4 $\degree C$ で遠心分離しました。mRNA 沈殿物を 70 % エタノールで 2 回洗浄して空気中で乾燥させてから、nuclease-free water に溶解させ、Agilent 2100 バイオアナライザ機器と Agilent RNA 6000 ナノキット (p/n 5064-1511) を使用して定量しました。次に、mRNA を 5 µg の量だけ作成して $-80 \degree C$ で冷凍しました。

次の2つの mRNA キャッピングメソッドを使用しました。

- 1) ARCA による共転写キャッピング
- 2) ワクシニアキャッピング酵素による酵素キャッピング

ARCA キャッピングでは、4:1 の割合の ARCA:GTP を含む NTP 混合 物で mRNA サンプルを転写しました。酵素によるキャッピングでは、 Vaccinia Capping System キット (New England Biolabs M2080) を使用して 5 ~ 10 μ g の精製済み未キャッピング (キャップされてい ない) mRNA をキャッピングしてから、Monarch RNA クリーンアップ キット (New England Biolabs T2040) を使用して精製し、分析前に nuclease-free water に溶出させました。

部位特異的 RNase-H 切断

Lapham, J. et al.⁴ の説明に従って、溶液内 RNase-H 部位特異的切断 をしました。15 nt キメラ 2'-O-メチル RNA/DNA プローブを mRNA サ ンプルを相補するように設計し(図 1)、次の 2 つの条件下で RNase-H 切断を誘導するのに使用しました。(1) 1 時間の熱アニーリングプログラ ムの実施後、37 °C で RNase-H により 30 分間切断、または(2)50 °C で耐熱性 RNase-H により 30 分間切断。切断後、Agilent 2100 バイオ アナライザ機器と Agilent Small RNA キット(p/n 5067-1548)で視覚 化する前に、5['] キャップを含む長さ 40 nt または 50 nt のオリゴヌクレオ チドを遊離し、分解したサンプルを Monarch RNA クリーンアップキット (New England Biolabs T2040)を使用して精製しました。

5' キャッピングされたオリゴヌクレオチドの LC-DAD/MS 装置構成

- 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。毎日 の実験を開始する前に、LC システムとカラムに 50 % MeOH + 0.1 % ギ 酸溶液を 30 分間通液し、アルカリ金属付加物をさらに減少させました。

分解した mRNAサンプルを、AdvanceBio オリゴヌクレオチドカラム (2.1 × 50 mm、2.7 μ m、120 Å、p/n 659750-702) で分離しました。 移動相と LC グラジエントを表 1 に、質量分析計の設定を表 2 に示しま す。MassHunter BioConfirm 10.0 と表 3 に示すデコンボリューション 設定によりデータ解析を実施しました。

Agilent 1290 Infinity II LC システム			
カラム	AdvanceBio オリゴヌクレオチド、2.7 μm、2.1 × 50 mm、120 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	時間 (分) % B 0 10 4 30 25 51 25~28 90		
カラム温度	50 °C		
流量	0.4 mL/min		
注入量	20 µL		
注入ごとの mRNA	2 µg		

表1.移動相とLC グラジエントの条件

表 2. 質量分析計の設定

Agilent 6545XT AdvanceBio LC/Q-TOF				
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	25 psig			
シースガス温度	275 °C			
シースガス流量	10 L/min			
V _{cap}	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			
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

MassHunter BioConfirm 10.0 の設定			
抽出クロマトグラム (MS)	ダイオードアレイ検出器		
MS1 遅延時間	0.06 秒		
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	10,000 ~ 30,000 Da		
質量ステップ	0.5 Da		
限られた m/z 範囲の使用	1,040 ~ 3,200		

結果と考察

通常、標準の LC/MS 分析では完全長 mRNA は大きすぎるため、最初 に RNase-H 部位特異的切断を最適化しました。5' UTR 内の別の配列 を相補するように 2 つの RNA/DNA キメラプローブを設計し、RNase-H 切断⁴を実行してヌクレオチド 40 または 50を作成しました (図 1A)。こ れにより、切断の完了時に 5' キャッピング構造を保持している長さ 40 nt または 50 nt のオリゴヌクレオチドを遊離させました。これらの小さい 5' キャッピングされたオリゴヌクレオチドは LC/MS 分析に適しています。



図 1. mRNA サンプル前処理と分析。(A) mRNA サンプルを相補するキメラプローブは、40 nt (紫) と 50 nt (金) で RNase-H 切断を実行するように設計しました。 下線が付けられたプローブヌクレオチドは DNA で構成し、それ以外は 2'-O-メチル RNA で構成しました。(B) 非耐熱性(上側の手法)または耐熱性 RNase-H (下側の手法)を使用したサンプル前処理法。下側の手法を使用した場合の合計分析時間:切断(30分) + クリーンアップ(10分) + LC/MS(25分) + データ解析 (10分) = 75分。

別の文献⁶で報告されているように、このような実験の一般的なワーク フローでは、非耐熱性 RNase-H 酵素を加える前に別の 1 時間の熱ア ニーリングを行い、37 ℃ でサンプルを切断します(図 1B、上側の手 法)。熱アニーリング時に使用する温度が高温(~95 ℃)でない場合 は、次の別の手順が必要になります。(1) 非耐熱性 RNase-H を変性し、 (2) 反応バッファに存在する二価カチオンにより mRNA 加水分解を促 進します。⁶

GC 含量が十分に高いキメラプローブを使用しているという条件におい て、適度に高い温度 50 ℃ で耐熱性 RNase-H により切断反応を促進 できると仮定しました (図 1B、下側の手法)。原理的には、高温では mRNA 二次構造が解かれることが予想されるため、高 GC 含量のキメラ プローブを相補的な mRNA 領域に結合して、耐熱性 RNase-H で切断 が可能となります。 図 2 に示すように、2 つの異なる高 GC 含量プローブ (プローブ 40 とプ ローブ 50、それぞれが > 60 % GC) と非耐熱性または耐熱性 RNase-H を使用して、部位特異的切断を行いました。生成した分解物を Agilent 2100 バイオアナライザ自動電気泳動ツールと Agilent Small RNA キット (p/n 5067-1548) を使用して分離し、仮想ゲルイメージにより定性的 に評価しました。

プローブ 40 とプローブ 50 はともに、すべての条件下で mRNA サンプ ルを正常に切断しました。ただし、プローブ 50 は非耐熱性 RNase-H を 使用した際に標的以外と結合し、好ましくない物質が生成されました(図 2、赤の矢印)。プローブ 50 は、耐熱性 RNase-H を使用した際に特異 的な切断をし、短期間でよりクリーンなサンプルを取得できました。



図 2. 耐熱性および非耐熱性 RNase-H を使用した部位特異的切断の評価。 プローブ 40 とプローブ 50 により、 ヌクレオチド 40 と 50 それぞれにおいて酵素に よる切断をしています。 赤の矢印は、 標的以外を切断した生成物を表しています。 L: ラダー。

次に、プローブ 50 + 耐熱性 RNase-H を使用して分解した未キャッ ピング mRNA を IP-RP LC/MS で分析しました。図 3 に、Agilent AdvanceBio オリゴヌクレオチドカラム (2.1 × 50 mm、2.7 μm、 120 Å、p/n 659750-702) において、未キャッピング 5' オリゴヌクレ オチドをプローブ 50 と残りの mRNA から分離した結果を示します。 MassHunter BioConfirm 10.0 に組み込まれている自動データ解析 ワークフローを使用して、UV クロマトグラムで同定したピークから質 量スペクトルを抽出してから、表 3 に示す設定により自動的にデコンボ リュートしました。表 4 に、各反応に存在する可能性があると予想される 未キャッピング、キャッピング中間生成物、および完全にキャッピングされ た種を示します。

未キャッピング mRNA は不安定な 5' 末端三リン酸基で構成されている ため、反応条件によっては、いくらかの量の加水分解された二リン酸また は一リン酸種3 が観察される場合があります。図4に、三リン酸および二 リン酸のピークから抽出された質量スペクトルとデコンボリュートした質 量を示します。最適化された耐熱性 RNase-H 切断プロトコルにより生成されたのは、少量(4.2%)の二リン酸種のみであり、一リン酸種は検出されませんでした。これは、反応条件が適切に穏やかであったことを示しています。

未キャッピングのサンプルでは、約15%の三リン酸オリゴヌクレオチド が、追加の非テンプレート化+Gヌクレオチドを含む配列変異体として検 出されました(図4D)。これはおそらくT7転写ずれが原因で発生してお り、転写配列の最初に反復Gヌクレオチドが存在する場合に一般的に観 察されます。⁷図5に示すように、ずれ配列変異体は、Cap1への次のメ チル化反応の前後で、ARCAで共転写キャッピングされたmRNAサンプ ル中にも確かに存在していました。ずれ配列変異体は明確な分離ピーク として同定できましたが、ベースライン分離されませんでした。キャッピン グされたオリゴヌクレオチドはすべて、未キャッピングのオリゴヌクレオチ ドから適切に分離されました。

表 4. 未キャッピング、キャッピング中間生成物、およびキャッピングされた種。赤の文字は、「p」の記号で示されるリン酸を含む 5' 末端ヌクレオチドを表しており、これは同化されるか、または酵素反応によりキャッピング構造に置き換えられる場合があります。

未キャッピング		キャッピング中間生成物		キャッピングされた種	
名前	配列	名前	配列	名前	配列
三リン酸	pppGGGGCC…	ニリン酸	ppGGGGCC…	ARCA Cap 0	3'-O-Me-m7GpppGGGGCC····
		G-Cap	GpppGGGGCC…	ARCA Cap 1	3'-0-Me-m7GpppmGGGGCC···
				Cap 0	m7GpppGGGGCC···



図3.5' 未キャッピングオリゴヌクレオチドのキメラプローブとサンプルマトリックスからのクロマトグラフィーによる分離。 ピーク1:5' ニリン酸、ピーク2:5' 三リン酸オリゴヌクレオチド

図 4.5' 未キャッピング mRNAオリゴヌクレオチドの質量スペクトル。図3のピーク1の(A) 抽出された質量スペクトルと(B) デコンボリュートした質量スペクトル。 図3のピーク2の(C) 抽出された質量スペクトルと(D) デコンボリュートした質量スペクトル。赤の数字は、二リン酸および三リン酸オリゴヌクレオチドに 起因する抽出された質量スペクトル内の各ピークの荷電状態を表しています。アスタリスクが付けられた質量ピーク(16131.16 Da と 16211.79 Da)は、推定される ニリン酸および三リン酸同定(配列挿入)と質量エラー < 10 ppm で一致していました(表 5)。パネル D では、三リン酸 + G 配列変異体(+345.17 Da)が、 三リン酸未キャッピングオリゴヌクレオチドと共溶出していました。

図 5. キャッピングされたオリゴヌクレオチドの未キャッピングのオリゴヌクレオチドからの分離。ピーク1~3:ARCA Cap 0、ピーク4~6:ARCA Cap 1。 キャッピングされた種には、0~2の非テンプレート化Gヌクレオチドがずれ配列変異体として含まれています。

図 6 に、ARCA Cap 0 と ARCA Cap 1 ピークから抽出された質量スペ クトルとデコンボリュートした質量を示します。報告されている文献²と同 様に、ARCA Cap 0 共転写キャッピングの効率は高い(91.3 ±1.76 %) ことがわかりましたが、完全ではありませんでした。その後 ARCA Cap 1 にメチル化されたサンプルでは、キャッピング効率は ARCA Cap 0 とそ れほど異なってはいませんでしたが(90.6 ±1.53 %)、これは実質的に すべての Cap 0 構造が完全に Cap 1 に変換されたことを示しています。

最後に、LC/MS メソッドを使用して、ワクシニア酵素キャッピングの効率 を評価しました。この反応は ARCA 共転写キャッピングより効率的であ り、サンプル配列に比較的依存しないと述べられていましたが³、今回の 実験でそうではないことが証明されました。メーカー推奨の反応条件を 使用したワクシニア酵素キャッピングでは、未キャッピングの物質、キャッ ピング中間生成物、および完全にキャッピングされたサンプルの混合物 が生成されました(図7と表 6)。図8に示すように、大部分のサンプル (90.9 ±0.6 %)は5'末端ニリン酸または三リン酸では未キャッピング のままであり、6.5 ±0.7 % はキャッピングされましたがメチル化されま せんでした(つまり、G-Cap)。キャッピングされて Cap 0 にメチル化さ れたのは、サンプルのわずか 2.6 ±0.5 % でした。 表 5. ARCA でキャッピングされたオリゴヌクレオチドおよびキャッピング中間生成物 のオリゴヌクレオチドの質量精度

ARCA 共転写キャッピング			
同定	理論質量	観察値の質量	精度(ppm)
ARCA Cap 0	16504.59	16504.62	2.01
ARCA Cap 0 + G	16849.80	16849.69	6.37
ARCA Cap 0 + 2G	17195.01	17195.03	1.50
ARCA Cap 1	16518.59	16518.85	15.77
ARCA Cap 1 + G	16863.80	16864.11	18.47
ARCA Cap 1 + 2G	17209.01	17208.73	16.13
ニリン酸	16131.52	16131.37	9.21
三リン酸	16211.51	16211.50	0.38
三リン酸 + G	16556.55	-	-

表 6. 酵素によりキャッピングされたオリゴヌクレオチドおよびキャッピング中間生成物のオリゴヌクレオチドの質量精度

酵素によるキャッピング				
同定	理論質量	観察値の質量	精度(ppm)	
Cap O	16490.59	16490.65	3.55	
Cap 0 + G	16835.80	16836.00	11.97	
G-Cap	16476.57	16476.65	4.85	
ニリン酸	16131.52	16131.49	1.95	
三リン酸	16211.51	16211.61	6.24	
三リン酸 + G	16556.72	16556.97	15.33	

図 6. ARCA Cap 0 mRNA オリゴヌクレオチドの質量スペクトル。(A ~ C) 図 5 のピーク 1 ~ 3 のデコンボリュートした質量スペクトル。(D ~ F) 図 5 のピーク 4 ~ 6 の デコンボリュートした質量スペクトル。アスタリスクが付けられた質量ピーク(16504.62 Da、16849.69 Da、17195.36 Da、16518.85 Da、16864.11 Da、および 17208.73 Da) は、推定される同定(配列挿入)と質量エラー < 20 ppm で一致していました(表 5)。緑の下線が付けられた文字は、T7 転写のずれにより非テンプレート化 された可能性の高いヌクレオチドを表しています。

図7. ワクシニア酵素でキャッピングされたオリゴヌクレオチドとキャッピング中間生成物のLC/MSの結果。(A) ピーク1~2: Cap 0 および +G 配列変異体、 ピーク3:G キャッピングされたオリゴヌクレオチド、(B~D) キャッピングされたオリゴヌクレオチドのデコンボリュートした質量スペクトル。アスタリスクが付け られた質量ピーク(16,490.60、16,836.00、および16,476.65 Da)はすべて、推定される同定(配列挿入)と質量エラー < 12 ppm で一致していました(表 6)。 緑の下線が付けられた文字は、T7 転写のずれにより非テンプレート化された可能性の高いヌクレオチドを表しています。

図8. LC/MS により導かれたワクシニア酵素キャッピングの最適化の結果

キャッピング効率を向上させるために、インプット mRNA の量を 0.5x および 0.25x に減少させました。これにより mRNA と比較して SAM、 GTP、およびワクシニアキャッピング酵素のモル比が増大しています。今 回の手法は比較的成功しており、未キャッピングの物質が減少した(0.5x = 73.6 ±1.3 % が未キャッピング、0.25x = 58.6 ±0.8 % が未キャッピ ング) 一方で、キャッピングされた物質が増大しました(0.5x = 17.7 ± 1.8 % が Cap 0、0.25x = 32.9 ±0.4 % が Cap 0)。単離された個々の 反応物質の濃度を増大させても、キャッピング効率は向上しませんでした (データは示していません)。これは、個々の成分が不完全であったこと を示しています。

結論

今回の実験では、mRNA 5' キャッピングを定量するためのより高速な LC/MS メソッドを開発しました。このワークフローで主に強化した点の 1 つは、耐熱性 RNase-H と GC 含量が十分に高いキメラ RNA/DNA オ リゴヌクレオチドプローブを組み合わせて使用し、別の熱アニーリングの 工程を実施しなくても mRNA 切断を促進したことです。GC 含量が極端 に高いプローブは、標的以外と結合するかまたは二次構造を形成する可 能性が高くなるため、適していない場合があります。実際、37 ℃ で切断 をした場合、プローブ 50 (GC 含量 67 %、図 2) において標的以外の切 断が発生しました。標的以外との結合は低反応温度で起こりやすいため、 高温で切断すると反応が促進されると同時に、よりクリーンなサンプルが 生成されました。別の実験では、耐熱性 RNase-H 切断する際、低 GC 含量 (40 %)のプローブでは成功しませんでした(データは示していま せん)。

次の2とおりの異なるキャッピング手法を使用して、今回のワークフロー を実証しました。(1) ARCA による共転写キャッピング、(2) ワクシニア キャッピング酵素による酵素キャッピング。ARCA キャッピングの結果は 文献とよく一致していましたが、ワクシニア酵素キャッピングによる今回 の結果は期待と異なるものでした。サンプル配列の分析では次の原因が 考えられます。つまり、最初の12種類のヌクレオチドはGC含量が90 %を超えて非常に高く、ワクシニア酵素キャッピングと干渉する場合があ る二次構造を形成する可能性が高かったということです。ホモポリマーG 反復では、非テンプレート化Gヌクレオチドを配列異性体として含める 場合もあるため(図5)、これらの結果からわかることは、この配列の5' UTR を再設計することにより、処理をより簡素化して生成物の品質を向 上できるということです。

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 近年、M.Beverly ら⁵によって同様の分析方法が発表されており、ここで はキメラプローブを磁気ビーズに共有結合することによって、RNase-H と切断したオリゴヌクレオチドの親和性の強化による部位特異的同時 切断が可能になりました。本実験では、非常に多くの項目を最適化した にもかかわらず、サンプルの特異的切断をすることができませんでした (データは示していません)。その理由は、サンプルの 5'末端を相補する キメラプローブの GC 含量が非常に高かったため、広範囲で非特異的結 合を引き起こしたことでした。サンプル中の分析困難な領域を回避するよ うにキメラプローブを設計できるため、今回のワークフローは柔軟性の高 いものであることがわかりました。さらに、シリカベースのスピンカラムを 使用したサンプルクリーンアップは、親和性ベースのアプローチと比較し て配列に依存することが少ないと考えられるため、サンプル回収率が増 大することが考えられます。

結論として、このアプリケーションノートでは、mRNA 5' キャッピングの 高分解能の高速分析において Agilent 6545XT AdvanceBio LC/Q-TOF が有効であることを示しました。また、本ワークフローは、医薬品開発ラ ボと品質管理ラボの両方において有用であると示唆されます。

参考文献

- Hassett, K. J. *et al*.Optimization of Lipid Nanoparticles for Intramuscular Administration of mRNA Vaccines.*Mol.Thera. Nucleic Acids* 2019, *15*, 1–11.
- Stepinski, J. *et al.*Synthesis and Properties of mRNAs Containing the Novel "Anti-Reverse" Cap Analogues 7-methyl(3'-O-methyl)GpppG and 7-methyl(3'-deoxy)GpppG. *RNA 2001*, 7, 1486–1495.
- Fuchs, A. L. *et al*. A General Method for Rapid and Cost-Efficient Large-Scale Production of 5' Capped RNA.*RNA* 2016, 22, 1454–1466.
- Lapham, J. et al.Site-Specific Cleavage of Transcript RNA. Methods in Enzymology 2000, 317, 132–139.
- 5. Beverly, M. *et al*.Label-Free Analysis of mRNA Capping Efficiency using RNase H Probes and LC-MS.*Anal.Bioanal. Chem.***2016**, *408*, 5021–5030.
- AbouHaidar, M. G. *et al*.Non-Enzymatic RNA Hydrolysis Promoted by the Combined Catalytic Activity of Buffers and Magnesium Ions.*Z Naturforsch C J Biosci* 1999, *54*, 542–548.
- 7. T.Conrad *et al*.Maximizing Transcription of Nucleic Acids with Efficient T7 Promoters.*Nat. Comm.Biol.***2020**, *3*, 1–8.

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, September 21, 2021 5994-3984JAJP PR7000-3098

