アプリケーションノート

# qPCRアッセイの最少化と 検量線作成の簡素化

**キーワード:**Multidrop Pico 8 Digital Dispenser、qPCR、 RT-PCR、液体分注、最少化、QuantStudio、直接混合、検量 線、微量qPCR反応

#### 概要

この試験では、Thermo Scientific™ Multidrop™ Pico 8 Digital DispenserとPicoIT 8 ソフトウェアを使用して、qPCR アッセイが最少化できることを実証および実践しました。 PicoIT 8 ソフトウェアは、簡単な3つのステップで、検量線に必 要な試薬のタイトレーションを含むqPCRアッセイプレートの セットアップを実現します。次いで、Multidrop Pico 8 Digital Dispenserを用いて、384ウェルのqPCRアッセイプレート に微量の試薬 (反応容量 5 µL)を分注して混合し、Applied Biosystems™ QuantStudio™ リアルタイムPCRシステム で増幅を行う準備をします。試験結果は、Multidrop Pico 8 Digital DispenserとPicoIT 8 ソフトウェアを利用することで、 最少量のApplied Biosystems™ TaqMan® qPCRアッセイプ レートで簡単にセットアップでき、input cDNAの良好な増幅が 得られることを示しています。

#### はじめに

Multidrop Pico 8 Digital Dispenserは非接触型の液体ディスペンサーで、アッセイ最少化とピコリッターレベルの溶液の分注 作業を可能にします。Multidrop Pico 8 Digital Dispenserは、 CV% (変動係数) 8%以下という正確さでわずか 11 pLの溶液 を分注可能で、96、384、1,536ウェルフォーマットのマイクロプ レートに対応しています。非接触型の分注テクノロジーは、試薬 のコンタミネーションリスクを排除します。さらに、Multidrop Pico 8 Digital Dispenserは分注した試薬を直接ミキシングす ることで、均一な反応を保証することができます。



Multidrop Pico 8 Digital Dispenserに搭載されたPicoIT 8 ソフトウェアは、使いやすいユーザーインターフェースを実現し ています。PicolT 8 ソフトウェアには、qPCRレイアウト用とし てプリセットされたプロトコルテンプレートが含まれています。 また、PCR functionを利用して簡単に自在にプロトコルを作成 することもできます。 プリセットのプロトコル gPCRテンプレー トの場合、ソフトウェアは3つのステップで進みます:まずプリ セットテンプレートを開き、お客さまのアッセイに合わせてプロ トコルテンプレートをカスタマイズして、新しく作成されたプロ トコルをMultidrop Pico 8 Digital Dispenserで実行します。 さらに、ソフトウェアは適切なフォーマット (.DA.csv) で関連 レポートを自動的に作成します。このレポートは、QuantStudio リアルタイムPCRシステムにインポートできます。サンプル情報 を含んだプレートレイアウトを直接インポートすることによっ て、QuantStudioでのプレートのセットアップ再作成のステップ が不要となり、qPCRのRunをすぐに開始できます。



TaqMan<sup>®</sup> qPCRアッセイのセットアップにおいてよくあるハードルの1つは、少ない反応容量でピペッティング精度を保証することです。 qPCRアッセイ開発では多くの場合、ピペッティングの誤差を理由として、やや多めの20 μL程度に反応容量を設定しています。精度を維持しつつ qPCRの反応容量を減らすことができれば、費用対効果が向上し、処理能力も改善されます。

この試験では、Multidrop Pico 8 Digital DispenserとPicoIT 8 ソフトウェアを利用して、384ウェルプレートに反応容量 5 µLのqPCRアッセイをセットアップすることに主眼を置いてい ます。この試験の各テクニカルリプリケートから得られたCt値 のCV%は、微量でのTaqMan®反応系の作成におけるPico 8 Dispenserの精度に関するエビデンスとなります。サンプル情報を含んだプレートレイアウトは、サーマルサイクルステップが 実行されるQuantStudio 5に直接転送されます。この試験の目 的は、Multidrop Pico 8 Digital DispenserとPicolT 8 ソフト ウェアを用いることで、qPCRアッセイを最少化でき、良好な検 量線を作成できることの実証です。これはqPCRワークフロー の一例であり、参考資料としてご利用いただくことはできます が、サンプルの種類、試薬、条件は実験デザインによって異なる ことにご注意ください。

#### 材料

#### 表1:このqPCRワークフローの例で使用する機器、試薬、消耗品

-	
機器	カタログNo.
Multidrop Pico 8 Digital Dispenser	Thermo Scientific, 5840600
QuantStudio 5、384ウェル用ヒートブロック	Applied Biosystems, A28575
E1-ClipTip電動シングルチャンネルピペット	Thermo Scientific, 4670000BT, 4670020BT, 4670040BT
試薬・消耗品	
TaqMan Gene Expression Master Mix (2x)	Applied Biosystems, 4370048
RT-PCR grade, Nuclease-free water	Invitrogen, AM9935
Triton X100、1%	Invitrogen, HFH10
TaqMan Assay A (20x, custom) FAM-MGB	Applied Biosystems, custom
TaqMan Assay (20x) RNaseP, VIC-TAMRA	Applied Biosystems, custom
cDNA (34 ng/µL)	Generated from Total RNA Control (human), Applied Biosystems, 4307281
384-well MicroAmp EnduraPlate、バーコード付	Applied Biosystems, 4483285
MicroAmp Optical Adhesive Film	Applied Biosystems, 4360954
ClipTipフィルターチップ (各サイズ)	Thermo Scientific, 94420043, 94420318, 94420813
4 ウェル 200 µL 分注ヘッドカセット for Pico 8	Thermo Scientific, LTR0003
8 ウェル 20 µL 分注ヘッドカセット for Pico 8	Thermo Scientific, LTR0004

#### 方法

まず、PCR反応コンポーネントを調製しました。Thermo Scientific™ E1-ClipTip™ ピペットを用いて、TaqMan® Assay 試薬を調製し、ヌクレアーゼフリー水でcDNAテンプレートを 希釈しました。Multidrop Pico 8 Dispenserで分注しやすくす るため、両方の混合液に最終濃度が0.1%となるようにTriton X100を添加しました<sup>1</sup>。表2は、反応容量 5 µLのTaqMan<sup>®</sup> qPCRアッセイにおける各反応コンポーネントの容量を示しています。

#### 表2:各ウェル (総量5 μL) のPCR反応コンポーネント

TaqMan Gene Expression Master Mix (2x)	2.5 μL
TaqMan Assay (20x) + Triton X100 (1%)	0.275 μL (0.1% Triton X100)
cDNA Template+ Nuclease free water + Triton X100 (1%)	2.225 μL (0.1% Triton X100)
総量	5 μL

本qPCRアッセイプレートの準備に用いた3ステップ:

プログラム済みの プロトコルテンプ レートを選択する

プロトコルをカスタマイズして保存 する (検量線などのパラメーターを 調整する)

分注プロトコルを実行し、 ソフトウェアのプロンプトに従う qPCRプレート の準備完了



#### ステップ1:

PicolT 8 ソフトウェアでプロトコルテンプレートを開きます。 File > Open Template > TaqMan Gene Expression MMx 384-wellの順にクリックします(図1、2)。



**図1. Multidrop PicolT 8 ソフトウェアのスクリーンショット** TaqMan Gene Expression Master Mix、384ウェル反応容量 5 µL、Multiplex のプロトコルテンプレートの選択肢を赤く囲んでいます。

#### 表3. タイトレーションの各レベルでPico 8 によって分注される容量

レベル	cDNA量
1	1.9 µL
2	0.95 µL
3	0.475 μL
4	0.238 µL
5	0.119 µL
6	59.4 nL
7	29.7 nL
8	14.8 nL
9	7.42 nL
10	3.71 nL
11	1.86 nL
12	0.928 nL
13	464 pL
14	232 pL
15	116 pL

#### ステップ2:

PicoITソフトウェアのPCR functionでプロトコルテンプレート をカスタマイズします。まず、リボンメニューからPCR function (図2の赤枠)を選択し、PCR functionにしたがって進めます。 Plate > Fluids > Unknown Groups > Standards/NTCs > Layout > Summaryの順にクリックして、実験デザインに固有 のパラメーターを設定します (図3、4、5)。サンプル数やリプリ ケートなどのカスタマイズも手動で行うことができます。この 試験では、PCR functionを用いて、NTCを4として一連のタイ トレーション曲線 (測定15ポイント、cDNA量 1.9 µLから始ま る1:2の線形タイトレーション、4リプリケート)を作成しました (表3)。すべての反応は 5 µLにノーマライズされています。



図2. TaqMan Gene Expression MMx (384ウェル) のプロトコルテンプレートを表示しているMultidrop PicolT 8 ソフトウェアのスクリーン ショット 赤枠はPCR functionのアイコンを示します。

CR					
late Fluids Unknown Group	ps Standards / NTCs Layou	it Summary			
Specify Maste	r Mix and one or more Unknow	n groups. All Primers v	vill be combined with each San	nple in a group.	
1 Master Mix	Unknown Group 1		ษตัว Unknown Gi	roup 2 🗙	
TQMN Gene Expression MMx	Replicates per sample 1		Replicates per sample 1		
2.5 µL	Brimore	Complex	Drimore	Complex	
1,1E-05 µL minimum	_ Phineis	Samples	Primers	samples	-
	✓ TQMN Assay 1 - FAM	CDNA 1	TQMN Assay 1 - FAM	CDNA 1	- +
	TQMN Assay 2 - VIC	1 Sample	✓ TQMN Assay 2 - VIC	1 Sample	
	1 Primer	0.5 µL	1 Primer	0.5 µL	
	0.275 µL	1.1E-05 µL minimum	0.275 µL	1,1E-05 µL minimum	
	1.1E-05 µL minimum		1.1E-05 µL minimum		
				Paula	

図3. PCR functionウィンドウと、Unknown Groupsタブにおけるこの試験のパラメーターセットのスクリーンショット

		Specify optional Standards and NTCs.	
✓ SI	tandards 🗸 St	andards/NTCs on each plate	✓ NTCs
Each Primer will be combined s Primers Standards ✓ TQMN Assay 1 - FAM ✓ TQMN Assay 2 - VIC 1 Standard 2 Primers 0.275 µL 1.1E-05 µL minimum	Replicates 4   Titration levels 15   Specify titration using   Highest and lowest volumes   Highest volume   Lowest volume   Highest volume	Replicates per primer 4 Primers TQMN Assay 1 - FAM TQMN Assay 2 - VIC 2 Primers 0.275 µL 1.1E-05 µL minimum	
Garrel		Distribution 1:2 (50%) V 18 levels without exceeding limits Minimum value is 1.1E-05 µL	Back Next

図4. PCR functionウィンドウと、Standards/NTCsタブにおけるこの試験のパラメーターセットのスクリーンショット

PCR			
Plate Fluids Unknown Groups	Standards / NTCs Layout Summary		
	130 Wells 1 Plate		
	NORMALIZE NORMALIZE	Normalize wells	
	Normalize to highest total total volume. Normalize to specified total volume.	Randomize each plate	
	Normalize using fluid class Aqueous+Triton X100 V		
	Normalize Do not normalize		
	Press Finish to create the PCR plates.		_
Cancel		Back Finis	sh

図5. PCR functionのSummaryウィンドウと、この試験のノーマライズのスクリーンショット

#### ステップ3:

Runをクリックして、分注プロトコルを開始します。E1-ClipTipを用いて試薬とサンプルをセットし、Multidrop Pico 8 Digital Dispenserで384ウェルマイクロプレートに分注します(図6の赤枠)。各ウェルに分注される具体的な容量については表2を参

照してください。Runの完了後、PicoIT 8 ソフトウェアは自動 的にレポートを作成します (.DA.csvファイルを含む)。次いでこ のファイルはqPCR のRunのためにQuantStudio Design and Analysisソフトウェアv1.5.1にインポートされます。



図6. この試験に関するカスタマイズ後のプロトコルとパラメーター設定のスクリーンショット 赤枠はRunコマンドを示します。

Multidrop Pico 8 Digital DispenserとPicoIT 8 ソフトウェア を用いてqPCRアッセイプレートを準備した後、このプレートに ついてQuantStudio 5システムの標準モードでqPCRを実行し ました。.DA.csvファイルをインポートするには、QuantStudio Design and Analysisソフトウェアv1.5.1において、File > Import Plate Setupの順にクリックします(図7)。



図7. QuantStudio Design and Analysisソフトウェアv1.5.1におけるプレートのセットアップと、この試験のパラメーター設定のスクリーンショット

# thermo scientific

## 結果

Multidrop Pico 8 Digital DispenserとPicoIT 8 ソフトウェア を用いることで、cDNAの増幅からqPCRアッセイプレートの 良好な検量線を設定することができました(図8)。Runレポー トは、Pico 8が各タイトレーションレベルでcDNAを分注して いることを示しています(表3)。QuantStudio Design and Analysisソフトウェアによる解析データは、検量線(VIC)の傾き は-3.325、プライマー効率は99.9%であることを示しています。 具体的には、R2値が0.996であり、試薬の分注が精密に行われ たことを示しています。各希釈点におけるcDNAの4リプリケート から得られたCt値のCV%は0.11~0.61%、すべての点の平均Ct のCV%は0.27%であり、極めて高い精度を示しています。



図8. この試験の増幅プロットと検量線を表示しているQuantStudio Design and Analysisソフトウェアのスクリーンショット

#### 結論

この試験は、Multidrop Pico 8 Digital DispenserとPicolT 8 ソフトウェアが、最少のqPCRアッセイの準備に優れたソリュー ションであることを実証しています。PicolT 8にあらかじめプ ログラムされているプロトコルテンプレートとPCR function を使用すれば、簡単な3ステップでqPCRアッセイプレートを セットアップできます。この結果は、Multidrop Pico 8 Digital Dispenserを使用することにより、アッセイ結果を損なうこと なくTaqMan<sup>®</sup> qPCRアッセイ反応の総反応容量を 5 µLまで 最少化できることを示しています。Multidrop Pico 8 Digital DispenserとPicolT 8 ソフトウェアは、検量線に必要な試薬の タイトレーションも実施し、段階希釈に伴う手動ピペッティング を不要とします。このタイトレーションは、相対定量PCRアッセ イや遺伝子発現アッセイ、プライマー効率バリデーション、PCR 効率バリデーション等における検量線作成プロセスのスピード アップに役立ちます。 さらに、各希釈液は、非接触型分注テクノロジーによって個々のウェル内に直接調製されます。この機能により、手動ピペッ ティングと段階希釈につきもののコンタミネーションや技術的 エラーのリスクが低下します。

Multidrop Pico 8 Digital DispenserとPicolT 8 ソフトウェ アを同時に使用することで、アッセイ開発における生産性向上 が実現可能です。これらは手動ピペッティングを削減し、試薬使 用量を減らし、高精度の結果をもたらす高速かつ最少化された qPCRアッセイを容易にします。

#### 参照文献

1. Weyent RS, Edmonds P, Swaminathan B. (1990) Effect of ionic and nonionic detergents on the *Taq* polymerase. Biotechniques Volume 9. Issue 3. Pages 308-309.

# 詳細はこちらをご覧ください thermofisher.com/multidroppico

These products are intended for General Laboratory Use. It is the customer's responsibility to ensure that the performance of the product is suitable for customers' specific use or application.

研究用にのみ使用できます。診断用には使用いただけません。 © 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. 実際の価格は、弊社販売代理店までお問い合わせください。 価格、製品の仕様、外観、記載内容は予告なしに変更する場合がありますのであらかじめご了承ください。 標準販売条件はこちらをご覧ください。thermofisher.com/jp-to LHC022-A2012OB

@ThermoFisherJP

## サーモフィッシャーサイエンティフィック株式会社

(i) TEL:0120-753-670 ≥ info.LPG.jp@thermofisher.com
オーダーサポート TEL:03-6832-9260 FAX:03-6832-9261
営業部 TEL:03-6832-9270 FAX:03-6832-9271



f facebook.com/ThermoFisherJapan