# Thermal Cycler Dice Real Time Systemシリーズ

新型コロナウイルス検査のための操作マニュアル

## -SARS-CoV-2 Direct Detection RT-qPCR Core Kit (製品コード RC330A) Primer/Probe N501Y (SARS-CoV-2) (製品コード RC344A)専用-

このマニュアルでは、SARS-CoV-2 Direct Detection RT-qPCR Core Kit (製品コード RC330A)、Primer/Probe N501Y (SARS-CoV-2) (製品コード RC344A)を用いてリ アルタイム PCR を実施する際の操作方法を説明します。実験操作に関しては、本キットの 取扱説明書に従ってください。

また、本装置は研究用機器であり、医薬品医療機器等法に定められる医療機器ではありません。

※本製品を弊社リアルタイム PCR 装置 Thermal Cycler Dice Real Time System シリーズで ご使用になる場合には、装置にデフォルトで設定されている正規化補正を解除したのち 解析を行ってください。正規化補正を設定している場合と解除した場合では、増幅曲線の 形状や Ct 値にわずかに差が生じることがあります。

解除方法は巻末の「Appendix: Thermal Cycler Dice Real Time System シリーズの正規化 補正解除方法」をご確認ください。

<装置とソフトウェアの起動>

- 1 Thermal Cycler Dice Real Time System 本体の電源を ON にする。
- 2 コンピューターの電源を ON にする。
- 3 食品環境検査用ソフトウェアを起動する。

<ランファイルの作成とランの開始>

- 1 ランファイルを新規作成する。
  - 1.1 解析タイプから絶対定量を選択する。
  - 1.2 多波長検出にチェック√を入れる
  - 1.3 OK ボタンをクリックする。

新規測定	
277±C b / →	
時年4月少1 ノ	
測定者名	<測定者の選択> ▼ 編集
	<b>OK</b> キャンセル

- 2 反応条件設定画面で PCR 条件を設定する。
  - 2.1 検出フィルターの「FAM」と「Cy5」にチェック√を入れる (ROX のチェック√は外す)
  - 2.2 Hold のパターンを1つ追加する。

●検出フィルター ▼ FAM ■ ROX	Cy5 Spee	ed Fast	ssociation time 2.0	↓ sec	
パターン セグメント	Hold 1	2 Step 1	PCR 2		
100 - - - - 50 - - - - - - - - - - - - - - - - - - -					
サイクル数	1	4	.0		
温度 <mark>(℃)</mark>	95.0	95.0	60.0		
時間(分、秒)	00:30	00:05	00:30		
データ取得					
4		194-2	RT Hold 2 Step PCR 3 Step PCR 融解曲線分析 力スタム	<ul> <li>パターン注意加</li> </ul>	セグメント追加

- 2.3 1つ目の Hold は、52℃、5分の設定にする。
- 2.4 2つ目の Hold は、95℃、10 秒の設定にする。
- 2.5 2 Step PCR の条件が、95℃、5 秒と 60℃、30 秒である事を確認する。
- 2.6 2 Step PCR のサイクル数は 45 にする。
- 2.7 Speed の設定は、Fast を選択する。

絶対定量 Multiplex	検出フィルター ▼ FAM	Cy5 Spee	ed Fast 🔻 Dis	ssociation time 2.0	sec 🖈
サンプル設定					
反応条件設定	パターン	Hold	Hold	2 Ster	PCR
12/07/1182/2	セグメント	1	1	1	2
結果/解析	100 -				
	- - - 50 - - - - - - - - - - - - - - - - - - -				
	サイクル数	1	1	4	5
	温度 <mark>(℃)</mark>	52.0	95.0	95.0	60.0
	時間(分、秒)	05:00	00:10	00:05	00:30
	データ取得				<b>V</b>

※上図は、Thermal Cycler Dice Real Time PCR System III の設定例です。

■他のランファイルからの設定読み込み

以前と同じ PCR 条件でランを行う場合には、他のランファイルから設定を読み込む ことができます。画面右上の"反応条件読込み"ボタンをクリックすると、ランファイ ルを選択するブラウザが開きますので、目的のファイルを選択して"開く"をクリック します。PCR 条件の他に蛍光フィルターの選択 ("データ取得") なども読み込まれます。



 サンプル設定画面でサンプル情報を入力する (ラン途中またはラン終了後に行っても良い)。

										表示切替 ③ マーク	② 名利	ウェル情報 入力 補	
	1	2	3	4	5	6	7	8	9	10	11	12	
A													
в													
с													
D													
E													

3.1 画面右上の入力ボタンをクリックする。

- 3.2 該当するウェルを選択し、サンプルタイプを選択する。
  - NTC: 陰性コントロールSTD: 陽性コントロール
  - **UNKN**: 検査対象サンプル

絶対定量 Multiplex	検	出フィルター EAM	ROX	v5					
サンブル設定									
反応条件設定 結果 /解析	A	NTC Cy5	Cy5	Cy5	Cy5	Cy5			
	в	STD Cy5 0.00E+000	Cy5	ウェル情報 サンブル名		×			
	с	UNKN Cy5	Cy5	- サンブルタ ⊂ターゲット マーカ	KN 】複数 連続設定				
	D	UNKN Cy5	Cy5	レプリケーマーク	·N設定	連続設定			
	E	UNKN Cy5	Cy5	検量線計 0.00	連続設定				
	F		Cy5	0mi	」カスタ t	L 0.0			
		HNKN	1						

3.3 必要に応じてサンプル名を入力する(省略可能)。 表示切替の「名称」を選択すると次のような表示になる。

											表示切替 © マーク	<ul> <li>名称</li> </ul>	
	1	2	3	4	5	6	7	8	9	10	11	12	
A	NTC		ウェル情報	ウェル情報設定									
в	STD		サンブル4 サンブル5 - ターゲッ	ら いけ UN ◆設定 〔	KN 】複数	•							
с	UNKN 検体1		Dye [	Cy5 ▼									
D	UNKN 検体2		マーク ( - 検量線)	<none td="" ▼<=""><td>連続設定</td><td></td><td></td><td></td><td></td><td></td><td></td><td></td></none>	連続設定								
E	UNKN 検体3 UNKN		10x	<u>* 555</u> カス! * カス!	94 <b>0.0</b>								

■他のランファイルからの設定読み込み

以前と同じ条件でサンプル設定をしたい場合は、他のランファイルから設定を読み込 むことができます。画面右上の"読込み"ボタンをクリックすると、ランファイルを選択 するブラウザが開きますので、目的のファイルを選択して"開く"をクリックします。

表示切	替	ウェル情報		
○ マ-	ク 💿 名称	入力	補助	読込み

★テンプレートファイルの利用も可能です。

上記の PCR 反応条件設定、サンプル設定を行った状態のファイルを

「テンプレートファイル」としてデスクトップに保存しておくと便利です。



新規ランファイルを作成する際は、まずテンプレートファイルを開き、ファイルメニュ ーから「別名で保存」を選択し、適切な保存先とファイル名を入力して保存して下さい。 必要に応じて設定を変更したうえでランを開始します。

Ø	ファ	·イル(F) 編集(E)	セクション(S) 解	析(A)	機器(I) ユーザ-	-(U) 表示(W)	リモート(R) ヘル	プ(H)
D		新規(N)	Ctrl+N	8				
		開<(O)	Ctrl+0					
-		閉じる(C)				d East T Di	esociation time 20	<u>^</u> 800
		データ管理(M)	Ctrl+M					
	_	(星友(C)	Ctrl+C					
	ſ	別名で保存(A)	Ctrl+A		Hold	Hold	2 Step	PCR
	L	==:1 7./T)			1	1	1	2
		記込み(1)		100-				
		フルレポート作成	(R)					$\land$
		印刷(P)	Ctrl+P	-		/		
		終了(X)		-		/		
	_			50 -				
				-				
				_				
				0-				
			サイクル数		1	1	4	5
			温度(℃)		52.0	95.0	95.0	60.0
			時間(分、秒	)	05:00	00:10	00:05	00:30
			データ取得					

- ※テンプレートファイルは、使用する Thermal Cycler Dice Real Time System を 制御する PC のみで利用可能です。別の装置制御用の PC へのファイル移動は 避けてください。
- 4 反応条件設定画面でランを開始する。
  - 4.1 反応用のチューブ(またはプレート)を本体にセットする。
  - 4.2 画面右下の反応開始ボタンをクリックしてランを開始する。



<結果の解析>

解析パラメーターの確認

- 1 増幅曲線を表示させる(N501Y Mutant 検出系(Cy5))
  - **1.1** 検出フィルターの Cy5 ボタンをクリックする。
  - 1.2 データ解析から増幅曲線を選択する。
  - 1.3 表示セレクトで解析対象のウェルを選択する。



2 Positive Control (N501Y Mutant、501N Wild) · Negative Control の確認 (Cy5)



Positive Control (N501Y Mutant)

Positive Control (501N Wild)



#### Negative Control



## 2.1 データ解析からテキストレポートを選択する。

検出フィルター	FAM HE	X ROX Oy5										2画面	全画面
ウェル	▲ サンプルタイプ	レプリケートマーク検出フィルタ	Ct值(CP) Ctd	(SDM)	Cutoff(Ct)	標準サンプル源(定量値(CP)	定量値(SDM)	Tm #1	Cutoff(Tm)		:;	テキストレポート	V
A7	NTC	Cy5	-	-					-				_
G7	STD	Cy5	-	-		0.000E+000 -			-	3	标形式	ウェル	~
G8	STD	Cy5	29.45	29.79		0.000E+000 -				1	辰示項目	□ 解析条件 ☑ CP法データ	
										5	翻項目	☑ SDM法データ	_
											マーリンフル名 マーサンフル名 マーレフリター	ら イブ トマーク	Î
											□ レブリケー □ 検出フィリ □ 最終蛍光 □ 最終蛍光 □ 兄信(CP) □ 平均C(値	ト名 リター (値 (Raw) (値 (CP)	Ŷ
												表示セレクト	V

2.2 以下を確認する。

	Cy5
	(N501Y Mutant 検出系)
Negative Control	不検出
501N Wild Template	不検出
N501Y Mutant Template	Ct≦30

- 3 増幅曲線を表示させる(501N Wild 検出系(FAM))
  - 3.1 検出フィルターの FAM ボタンをクリックする。
  - 3.2 データ解析から増幅曲線を選択する。
  - 3.3 表示セレクトで解析対象のウェルを選択する。



4 Positive Control (N501Y Mutant、501N Wild)・Negative Control の確認 (FAM)



#### Positive Control (N501Y Mutant)

## Positive Control (501N Wild)







- 検出フィルター FAM HEX ROX Cy5 2画面 全面面 (SDM) Cutoff(Ct) 」サンプルタイプ ターゲットマーク レプリケートマーク 検出フィル 標準サンプル湯|定量値(CP) 定量値(SDM) Tm #1 Ct値(CP) Cutoff(T デーク解析 Ċιλ Ct テキストレポート NTC FAM ... ... ... ... ... ... G3 STD FAM 27.72 28.69 0.000E+000 表示形式 Ċιル G4 STD FAM 0.000E+000 ... ... ... ... □ 解析条件
   ✓ CP法データ
   ✓ SDM法データ 表示項目 詳細項目
- 4.1 データ解析からテキストレポートを選択する。

4.2 以下を確認する。

	FAM
	(501N Wild 検出系)
Negative Control	不検出
501N Wild Template	Ct≦30
N501Y Mutant Template	不検出

- 5 検体の結果確認
  - 5.1 検体の増幅を確認する。



- 5.2 データ解析からテキストレポートを選択する。
- 5.3 Ct 値(CP)を確認する。

検出フィ ター	FAM	X ROX	Cy5											2画面	全画面
ウェル	▲ サンプルタイプ	ターゲットマーク	レプリケートマー	検出フィルター	Ct値(CP) Ct	(SDM)	Cutoff(Ct)	標準サンプル濃 定	量値(CP)	定量値(SDM	) Tm #1	Cutoff(T	#! Z##	テキストレポート	~
B3	UNKN			Cy5									2 2000	1 44104	
B3	UNKN			FAM				-			-	-	表示形式	VILV	
C8	UNKN			Cy5	33.36	35.27		-		-	-	-			
C8	UNKN			FAM	-	-		-	-	-	-	-	表示項目	□ 時竹奈仟	
				_									Nin/TO	☑ SDM法データ	,
													<ul> <li>マウエル</li> <li>マウエル</li> <li>サンブルタ</li> <li>サンブルターゲット</li> <li>マレブリケー</li> <li>マレブリケー</li> <li>マ枝出27(1)</li> <li>最終量光</li> </ul>	5 イブ マーク A-マーク トマーク 小名 ター 値 信	*
<												>		表示セレクト	

#### 6 結果判定

以下のような判定を行う。

FAM (501N Wild 検出系)	Cy5 (N501Y Mutant 検出系)	判定
<b>Ct&gt;40</b> または不検出	Ct≦40	<b>N501Y</b> 変異あり
Ct < 10	Ct \ 10 またけて 按山	<b>N501Y</b> 変異なし
01≧40	01/40 または小俠山	(Wild Type(501N))
Ct>40 または不検出	Ct>40 または不検出	判定不能

 501N Wild 検出系、N501Y Mutant 検出系共に Ct>40 または不検出であった場合 は、判定不能です。予め PCR 検査で SARS-CoV-2 陽性であることを確認したサン プルを使用した場合、検出限界以下あるいは他の変異の可能性があります。SARS-CoV-2 の有無が未確認のサンプルを使用した場合、SARS-CoV-2 陰性の可能性が あります。 7 解析結果の出力

(増幅曲線)

1 出力したいサンプルを選択し、画面上で、右クリックする。

2 データ出力から Excel か、レポート出力から Word · PowerPoint を選択する。

検出フィルター FAM F		2画面 全画面
100		デーダ解析 増幅曲線 🔹
80-	24-	グラフ表示 Primary Curve ▼
	Xm +	解析設定
- <u>-</u>	Yea >	フィルター Cy5 ▼ 道用
40 -		ベースライン 閾値
20- 第 20- 第		Auto     Fluorescence Cycle     O    Auto
<sup>20</sup> 0 -	● 原出力 → データ出力 → CSV	Manual 9.84 ▼ 8 ▼
1 2 3 4 5	レポート出力 ・ Excel 21 21 22 23 24 25 26 27 28 29 30 31 32 33 34 35 36 37 38 39 40 41 42 印刷	43 44 45

(テキストレポート)

- 1 出力したいサンプルを選択し、画面上で、右クリックする。
- 2 データ出力から Excel を選択する。

ウェル	▲ サンプルタイプ	レプリケートマーク	検出フィルター	Ct値(CP)	Ct値(SDM)	結果(F)	判定(F)	結果(CP)	判定(CP)	結果(SDI	デーク解析	テキストレポート ▼
D5	UNKN		Cy5			. –	Nega.	-	Nega.	-	2 244	2.1231.020-1
D6	PC		Cy5	25.42	25.37	+	OK	+	OK	+	表示形式	<b>ウェル</b> ▼
D7	NC		Cy5				OK	-	OK	-		277+1 - + 2 /4
											表示項目	<ul> <li>         ・一 解析条件         ・・・・・・・・・・・・・・・・・・・・・・・・・・・・・</li></ul>
			32-								詳細項目	V SDM法データ
			ウェルの並べ方	•							✓ ウェル ■ サンプルタ	
			31/2/10/1E0/12	<b>庆</b> 9							▼ サンブルタ	17 🗐
			データ出力	•	CSV						▼ レブリケート	トマーク
					Excel						■ レノリク = I ▼ 検出フィル	~ー ター
											■ 最終蛍光	値(Raw)
											■ 最終国光	10 -
•									_	•		+ Ы
								_				表示セレクト 🔽

<ソフトウェアと装置の終了>

- 1 食品環境検査用ソフトウェアを終了させる。
- 2 コンピューターを終了させて、電源を切る。
- 3 Thermal Cycler Dice Real Time System 本体の電源を切る。

Appendix1: Thermal Cycler Dice Real Time System シリーズの正規化補正解除方法

- A. Thermal Cycler Dice Real Time System III の場合 (Software Ver. 3.01C/3.01D)
  - 1. ソフトウェア画面左上のユーザー(U) → 設定(S)をクリックする。



2. ユーザー設定内の解析タブを選択し、正規化補正のチェックを外す(赤矢印)。

Software Ver. 3.01C



3. 右下の OK をクリックしウインドウを閉じる。

- B. Thermal Cycler Dice Real Time System II/Lite の場合(Software Ver. 2.11C)
  - 1. ソフトウェア画面左上のユーザー(U) → 登録(M)をクリックする。

📝 食品環境検査用ソフ	トウエア	
ファイル(F) 機器(I)	[ユーザー(U)] ヘルプ(H)	
D 🗲 🔳 🏉 🛍	変更(C)  ▶ Ō 設定(S)	?
	登録(M)	

 新規 → 適当なユーザー名を入力(例:正規化補正 OFF) → 追加の順に操作し、 上部リストにユーザー名が追加されたのを確認したのち、下部の OK をクリックし ウインドウを閉じる。

登録		
名前	XE	
正規化補正C		
ユーザー	#T055	tr ta
	#IEOFF	新規
		Ť BR
	OK キャンセル	]

3. 新規 Run file を作成する際に、前項で登録したユーザー名を選択し OK をクリック。

新規測定		X
解析タイプ	絕対定量	▼ 🔽 多波長検出
測定者名	正規化補正OFF	▼ 編集
	<u>ОК</u> <i>キャンセル</i>	,

📝 食品環境検査用ソフトウエア - [NewDocument\_6] D 🗲 🖬 🎒 🗈 🗠 က ଲ 🔍 🔍 🕭 💡 変更(C) ۲ 設定(S) 検出フィルターー 絶対定量 Multiplex 登録(M) FAM HEX サンプル設定 2 4 1 3 5 6 反応条件設定 FAM FAM FAM FAM FAM FAM А 結果/解析

.....

4. ソフトウェア画面上部のユーザー(U) → 設定(S)をクリックする。

5. ユーザー設定内の解析タブを選択し、正規化補正のチェックを外す(赤矢印)。

|\_...

L....

ユーザ設定		
解析レポート作成	/印刷   グラフのプロパティ   テキストレポート   ランプ	全てリセット
解析 增幅曲線	☑ 各ターゲットごとにパラメータを設定	Utyk
	ベースライン 💿 Auto 💿 Manual	
	閾値 💿 Auto 🔘 Manual	N
▶ →	📃 正規化補正	
融解曲線	☑ 各ターゲットごとにパラメータを設定	
測定値分布	☑ 各ターゲットごとにパラメータを設定	
	閾値 💿 Auto 💿 Manual	
-		-
		N
۶Æ		OK         キャンセル

6. 右下の OK をクリックしウインドウを閉じる。これ以降は Run file を作成または解 析する際に、同じユーザー名を選択すれば、常に正規化補正が解除された状態とな る。

## C. 上記以外の Software Ver.の場合(III/II/Lite 共通)

**※Software Ver. 2.11C/3.01C/3.10A** 以外は Run file ごとに正規化補正を解除する 必要がある。

1. Run file を開いた状態で、解析(A)  $\rightarrow$  基本設定(S)をクリックする。



2. 正規化補正のチェックを外す(赤矢印)。

基本設定
ースムージング
Amplification Averaging Points 5
Dissociation Averaging Points 5
正規化補正
Amplification Plots 🔲 🖛
OK ++>>セル

3. 左下の OK をクリックしウインドウを閉じる。