CronoSTAR[™] 96 Real-Time PCR Systemシリーズ

新型コロナウイルス(SARS-CoV-2)検査のための操作マニュアル

-SARS-CoV-2 Direct Detection RT-qPCR Kit(RC300A, RC30JW)専用-

このマニュアルでは、SARS-CoV-2 Direct Detection RT-qPCR Kit (製品コードRC300A, RC30JW) 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。

CronoSTAR 96 リアルタイムPCR装置とソフトウェアの起動

- 1. リアルタイムPCR装置本体の電源をONにする。
- 2. コンピューターの電源をONにする。
- 3. CronoSTAR 96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. New Experimentをクリックし名前を入力してランファイルを作成する。



2. Run Settingタブの[+]を押し[Reverse Transcription],[2 Step Amplification],[Cooling]を追加する。



- 2.1. [Reverse Transcription]の設定
 - 2.1.1. Step1は、52℃、5分、Ramp 6.1℃/sの設定にする。
 - 2.1.2. Step2は、95°C、10秒、Ramp 6.1°C/sの設定にする。
- 2.2. [2 Step Amplification]の設定
 - 2.2.1. Step1は、95℃、5秒、Ramp 6.1℃/sの設定にする。
 - 2.2.2. Step2は、60°C、30秒、Ramp 6.1°C/sの設定にする。
 - 2.2.3. サイクル数を×40から×45に変更する。
- 2.3. [Cooling]の設定
 - 2.3.1 Step1は、35℃、30秒、Ramp 6.1℃/sの設定にする。



3. ExperimentのTube Type [Clear]、Reaction Volume [50µL]、Lid Heating[105℃]に設定する。



※ 他のランファイルからの設定を読み込む場合

User Name: user Switch User	New Experin	ment	X
Quick Start Recent Files		任意で変更可能	
>>New Experiment >>New Experiment From Existing Experiment	Experiment N	lame user 20210415083551	
>>Open Data File		∖変更不可	Connect
>>Instrument Management Default Instrument: CronoSTAR 96-6ch-#10		New	incei

以前と同じPCR条件でランを行う場合には、New Experiment From Existing Experimentを クリックし参照したいファイルを選択、名前を入力してランファイルを作成する。

- 4. Sample Setting タブでサンプルの情報を入力する(後で入力可)。
 - 該当するウェルを選択し、サンプル名、タイプを選択する。DyeはCy5にチェック√をいれる。

Negative: 陰性コントロール

Positive:陽性コントロール

Unknown: 検査対象サンプル



- 5. 装置にサンプルをセットする。
- 6. Run MonitoringタブのRunをクリックしてランを開始する。



<u>結果の解析</u>

1. Analysisタブで測定結果を確認する。



2.Analysis settingボタンを押すとウインドウが表示される。



3. Analysis setting ウインドウのNormalization Methodを選択する。OKをクリックして一旦閉じる。

Analytical Method: 🔘 Aut	o Threshold 🔘 Manual Thresho	old 💿 Normalization Method	Restore
Dye	Gene	Auto Threshold	Manual Threshold
Cy5		814.46	1195.45

4. Analysis setting ウインドウを再度開き、Baseline Gain Calibrationのチェックを外す。

Amplification F	Gene and Sample	
Analysis Mode:	🗌 Reference Dye 🔲 Baseline Gain Calibration 📗 Reverse Curve 🔲 Isothermal	
Baseline	All Selected Rows: Start Cycle: 🗧 茾 End Cycle: 📮	

5. Amplification Curve画面で増幅曲線を確認する。Cy5にチェック√が入っていることを確認。



自動判定では曲線が見られなくとも陽性と判断される場合があるので Raw Curve で必ず増幅曲線、曲線の立ち上がりについてコントロールと比較して確認する。

6. Result Table 画面の Ct 値を確認する。

Resul	t Statist	ics								
Well	Sample ID	Sample	Sample Type	Dye	Gene	Ct	Concentration	Concentration Unit	Standar	1
F8			Unknown	Cy5		25.454		IU/ml	-	*
G8			Unknown	Cy5		-		IU/ml	-	
H8			Unknown	Cy5		-		IU/ml	-	
A9			Unknown	Cy5		-		IU/ml	-	
B9			Unknown	Cy5		-		IU/ml	-	
C9			Unknown	Cy5		-		IU/ml	-	
D9			Unknown	Cy5		-		IU/ml	-	
E9			Unknown	Cy5		-		IU/ml	-	
-		1		~ ~		1			Þ	

結果の判定方法(詳細はキット(RC300A)の説明書を参照)

コントロール反応の判定

- ・陰性コントロールは、不検出であることを確認する。
- ・陽性コントロールは、Ct 値が30 以下であることを確認する。
 ※陰性コントロールでCt値が算出された場合には、「補足②」に従いCt値算出の原因を確認する。また、Ct値算出の原因としてコンタミネーションの疑いがある場合は、反応液の調製場所や器具類を除染したうえで再反応を行う。

検体の測定結果の判定

- ・Ct 値が40 以下の場合は、陽性と判定する。
- ・Ct 値が40 より大きい場合、または不検出の場合*は検出限界以下と判定する。
- ※検出限界以下と判定された場合でも、40 サイクルより後で増幅曲線の立ち上りが認められた 場合は、低コピーの新型コロナウイルス遺伝子が存在する可能性があります。

<u>解析結果の出力</u>

1. Result Table 画面の Export Excel をクリックして名前を付けて保存する。

							R	lesult Table		3
Result Statistics										
Well	Sample ID	Sample	Sample Type	Dye	Gene	Ct	Concentration	Concentration Unit	Standar	
F8			Unknown	Cy5		25.454	-	IU/ml	-	*
G8			Unknown	Cy5		-	-	IU/ml	-	
H8			Unknown	Cy5		-	-	IU/ml	-	
A9			Unknown	Cy5		-	-	IU/ml	-	
B9			Unknown	Cy5		-	-	IU/ml	-	
C9			Unknown	Cy5		-	-	IU/ml	-	
D9			Unknown	Cy5		-	-	IU/ml	-	
E9			Unknown	Cy5		-	-	IU/ml	-	*
				~ -					+	

<u>ソフトウェアと装置の終了</u>

- 1. 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. 本体画面上でシャットダウンを押す。
- 4. 本体の電源を切る。

【補足①】前処理(核酸の簡易抽出)をリアルタイム PCR 装置で実施する場合

- 1. New Experimentをクリックし名前を入力してランファイルを作成する。
- 2. Run Settingタブの[+]を押し[Initial Denaturation],[Cooling]を追加する。
- 3. Initial Denaturation Step1は、95°C、5分、Ramp 6.1°C/sの設定にする。
- 4. Cooling Step1は、4℃、30秒、Ramp 6.1℃/sの設定にする。



- 5. 装置にサンプルをセットする。
- 6. Start ボタンを押しランを開始する。

【補足②】判定の際の注意

本機器はRunが終了すると、自動解析によりベースライン補正を行います。サイクル初期は蛍光値が 揺らぎやすく、自動解析によりこの区間がベースライン範囲に指定されてしまうと、不適切な補正に より増幅していない反応があたかも増幅したような曲線に補正されCt値が算出される場合がありま す。

そのため、判定の際には、以下の手順にしたがって、上述の自動補正による影響の有無を必ずご確認 ください。



Amplification Curve の結果

サイクル初期で Ct 値が算出されている事例。Raw Curve でも増幅が見られるか確認する。





サイクル初期 10 サイクル目くらいまでは蛍光値が揺らぎやすいため、どの区間をベースライン範囲 として認識しているか確認する。 ベースラインの確認と変更方法



①Analysis Setting アイコンをクリックする。

②Analysis Setting の変更したい Well と Manual Baseline を選択する。

③Start Cycle と End Cycle の数値を変更して OK をクリックする。



ベースラインの変更後

End Cycle を4から10に変更するとCt値が算出されなくなる。