

Applied Biosystems 7900HT Fast Real - Time PCR System

Automation Controller 簡易操作ガイド

SDS software version 2.3





Rev.A

研究用にのみ使用できます。診断目的およびその手続き上での使用は出来ません。

The PCR process and 5' nuclease process are covered by patents owned by Roche Molecular Systems, Inc. and F. Hoffmann-La Roche Ltd.

Applied Biosystems is a registered trademark and AB (Design) and Applera are trademarks of Applera Corporation or its subsidiaries in the US and/or certain other countries.

詳細については、機器付属のユーザーガイド(英文)を必ず参照ください。

© 2006 Applied Biosystems Japan Ltd,. All rights reserved.

本ガイドはApplied Biosystems 7900HT Fast Real - Time PCR Systemの操作を行う際に必要となる部分のみを記載している簡易操作ガイドです。

詳細は装置付属の英文ユーザーガイドをご確認下さい。また本ガイドには知的財産に関するいかなる情報も含んでおりません。知的財産に関するする情報は英文ユーザーガイドに記載されておりますのでご参照ください。

SDSシステムソフトウェア簡易操作ガイド

Automation Controller使用方法 ~Zymark[®] Twister[®] Microplate Handler を使用した連続ラン~

<u>概要</u>

Applied Biosystems 7900HT Fast Real-Time PCR Systemの自動プレートハンドラ付属モデルでは Zymark[®] Twister[®]マイクロプレートハンドラーを使用した自動連続ランに対応することができます。 またAutomation Controller Softwareによりプレート情報をBatchにまとめて連続ランを行うことができます。

<u>連続ラン用Plateドキュメントの作成と連続ランの実行フロー</u>

自動プレートハンドラを用いて連続ランを行う前に、以下のいずれかの方法でプレートの情報を Automation Controller Softwareに登録(キューの登録)する必要があります。

- 1.プレート情報の登録
 - ・SDSソフトウェアを用いて作成した個別のファイルからPlate Queueを作成する方法 (p.4~) 数枚のプレートを使用するランで、プレート毎に設定が異なる実験の場合に利用します。
- ・ テンプレートファイルを使用してファイルを複製し、Queueに登録する方法 (p.6~)
 全プレートすべて同一の実験フォーマットでランを行う場合に利用します。
- 2. Automation Controller を用いた連続ランの開始 上記いずれかの方法で登録されたQueueの情報に基づいてランを実行します。
- 注: Automation Controllerを用いた連続ランを行う場合、バーコードつきプレートを必ずご使用ください。

個別のファイルからPlate Queueに登録する方法

- 1. SDS v2.3ソフトウェアでプレートドキュメントを作成します。
- ドキュメント作成についての詳しい情報は別冊の納品検収補足資料・簡易操作ガイドをご参照ください。
- 2. Tools メニューから Document Information を選択します。

0	SDS	2	2.3									
File	Edit	1	⊻iew	Tools	Instrument	<u>A</u> nalysis	<u>Window</u>	Help)			
		á	•	Det	ector Manage	er	Ctrl+T	2	>			
0	Aut	0	tes	Dve	ker Manager	12	Ctrl+J	20	tio	on		
	1	2	3	Tab	ole Settings M	lanager	Ctrl+K	4	15	16	17	18
A				Doc	cument Inform	nation	Ctrl+U					
в				Loc	al User Acco	ount Manag	er					
C		1		Cha	ange Passwo	ord						
Ě	-	-		Opt	ions		Ctrl+Q	F				
D	_	_								-	-	
E												

3. Document Information ダイアログが表示され、Barcode フィールド内にカーソルが現れます。

Plate Description	Information	Operator
File Name: Instrument: Assay: Last Modified: Acquisition started: Acquisition ended:	II Auto test Unknown Absolute Quantification 08/03/2006, 16:58:16	This name will appear on reports.
Data collected on: Plate Format:	Unknown 384 Wells Clear Plate	
The comments you e	nter here will appear on reports	
? Print		OK Cancel

4. 付属のバーコードリーダーを使用し、下図の要領でボタンを押してバーコードを認識させます。



5. Barcode フィールドに読み取られたバーコードが表示されます。Document Information 画面下部の OKボタンをクリックします。

Plate Description	
Barcode:	384N0185OR
File Name:	Auto test

6. Instrument タブから Queue を選択し、Send To Queue ボタンをクリックします。

rmal Cycler Real-Time Queue Send To Queue Collect vata Collection Stamp Pre:	Collect post-read data only
Send To Queue Collect	Collect post-read data only
Pre:	
Pre:	
Post:	
LOST.	

- 7. ドキュメントの保存確認のメッセージが表示されます。Yes ボタンをクリックします。
- 8. Queueに登録するメッセージが表示されますので、Yes ボタンをクリックします。 この操作により Plate Queue に情報が保存されます。
- 9. メッセージが表示されますので、OKを選択します。

10. 複数のプレートドキュメントを登録する必要があれば、上記1~8の操作を繰り返します。

11. すべてのプレートドキュメントを登録したら、ソフトウェアを終了します。

<u>テンプレートファイルからプレートドキュメントを複製し、Queueに登録する方法</u>

- 1. SDS v2.3ソフトウェアでプレートドキュメントを作成します。
- ドキュメント作成についての詳しい情報は別冊の納品検収補足資料・簡易操作ガイドをご参照ください。 2. 作成後、ドキュメントをテンプレートファイル(.sdt形式)で任意のフォルダに保存します。

5				×
C 2.3A	-	£	ď	
Auto 1.sdt				Save
SDS 7900HT Template Document (*.sdt)		-	0	Cancel
	S 2.3A Auto 1.sdt SDS 7900HT Template Document (*.sdt)	S Calculate Spectral Content (*.sdt)	S 2.3A C Auto 1.set SDS 7900HT Template Document (*.set)	S 2.3A C Auto 1.set SDS 7900HT Template Document (*.set) C C C C C C C C C

- 3. 2.から連続で操作する場合はすでにドキュメントが開いているので、次に進みます。 ドキュメントが開いていない場合は保存したテンプレートファイルをFile メニューからOpenで開きます。
- 4. Instrument タブから Queue を選択し、Send To Queue ボタンをクリックします。

iermai Cycler Real-Time Gueue	
Send To Queue	📕 Collect post-read data only
-Data Collection Stamp	
Pre:	
Post:	

5. Template Batchダイアログが表示されます。New ボタンをクリックします。

🗵 Templ	ate Batch			X
		Plate ID		
				=
1	D:\Applie	ed Biosystems\SDS	S Documents'(2.3Auto	
New	Open	Import		Clear
Plate Directory:	D: VApplied Biosyst	tems\SDS Documer	ntsVhashizume	
?	Drowse		Create	Done

6. バーコード入力ダイアログが表示されます。 付属のバーコードリーダーを使用し、下図の要領でボタンを押すとプレートのバーコードが読み取られ、 Template Batch ダイアログにバーコードが自動的に入力されます。この操作をプレート枚数分繰り返します。



手入力する場合はPlate IDにバーコードを入力してください。入力を繰り返すときは Create Another ボタンをクリックします。

🗵 Nev	/ Plates			×
Plate ID:	I			
		Create Another	ок	Cancel

- 7. すべてのプレートのバーコード読み取りが終了したら、New Plates ダイアログのOKボタンをクリックします。
- 8. 読み取ったバーコードがTemplate Batch ダイアログボックスに表示されていることを確認し、 Browse ボタンを押してプレートドキュメント保存先を指定します。

	Plate ID	
1 384N018	50R	<u> </u>
2 384N018	5PS	
3 304N010 4 384N018	581	
5 384N018	55V	
		T
New	Open	Clear
140.17	injort	0.04
Plate Directory:	D: Applied Biosystems/SDS Documents/2.3A	uto
e Directory:	D: \Applied Biosystems\SDS Documents\2.3A	uto

 Create ボタンを押すことで登録したバーコードのプレートドキュメントが複製され、 Queue に自動登録したことを示すダイアログが表示されます。OK ボタンをクリックします。



10. Template Batch ダイアログ下部のDone ボタンをクリックし、ダイアログを閉じます。 11. ソフトウェアを終了します。

<u>Automation Controller を使用した連続ラン</u>

Zymark[®] Twister[®] 7900HTシステム本体 Automation Controller





384 well プレート、Fast 96 well プレートおよび Optical capで蓋をした96 well プレートの場合 はそのままスタックにセットします。 Optical Adhesive Filmでシールした96 well プレートを使用する場合、MicroAmp® Snap-on Optical Film Compression Pad (PN 4333292)を右図のように組み合わせ、 スタックにセットします。

※ラン開始前に排出スタックにプレートが残されていないことをご確認ください。

<u>Automation Controller を使用した連続ラン</u>

1. SDS Automation Controller 2.3 ソフトウェアを起動します。



2. 7900HTシステム本体、バーコードリーダーおよびAutomation Controllerとのコミュニケーションを 自動的にチェックします。



- 3. すべてパスすると、Automation controller が起動します。画面構成は下図の通りです。
 - A:装置とのコミュニケーション、ランに関する情報が表示されます。
 - B: Queuelc登録されているプレートがリスト表示されます。未登録の場合は空欄になります。 Plate Queue Queuelc登録されているラン未実行のプレートIDが表示されます。 Plate Status 装置のブロック、サンプル温度変化を常にモニターしグラフ表示します。 Processed Plates ランが終了したプレートIDが表示されます。
 - C: ランの実行および使用するスタックに関する設定、自動解析設定の設定を行います。

SDS Automation Control	ler	
File Edit Instrument Tools Help		
2 4 4 4 7 2 8 8 4	· 🔍	
Instrument: Overview Status: Edite Block Type: 384 Wells Block Serial Number: 200027 Batch Status: Queued Plates: 0 Processed	Plate Queue Run Status Processed Plates 0 Plates Currently Queued Barcode Plate Name Barcode Barcode Bar	Run Type
	Save List Add Plates Yew Plate Remove Selected B Automation Robot Sattings Select the stads that contain the plates for this batch IF Stack 1 Stack 2 Stack 3 Stack 4 Restack plates after If Stack 2 Stack 3 Stack 4 Restack plates after If Stack 3 Stack 4 Restack plates after If Stack 3 Stack 4 Restack plates after If Stack 5 <	emove All
		tart Batch Idle 🧕

4. Automation Robot Settings内の使用するスタックにチェックを入れます。

E CDC Automation Controll		
SDS Automation Controll		
rie cut pistrulient 1008 Hep	a	
	4	
Instrument Overview	Plate Queue Run Status Processed Plates	
Status: Idle	0 Plates Currently Queued	
Block Type: 384 Wells Block	# Barcode Plate Name	Run Type
Serial Number: 200027		
Batch Status		
Queued Plates: 0		
Processed Plates: 0		
Batch Long Show Datale		
Silow Decais		
Currently Running Plate		
Run Time Remaining:		
Name:		
Run Type:		
	Save List Add Plates Yew Flate Remove Selected R	temove All
	Automation Robot Settings	
	Select the stacks that contain the plates for this batch	
	🗹 Stack 1 🔲 Stack 2 🔲 Stack 3 🗐 Stack 4 📄 Restack plates after run	
	Enable Batch Settings	
	T Auto-Analysis Workflow T E-mail notifications	ettings
	Open/Close Door	Start Batch
		Idle 🧝

※オプション

Enable Batch Settings では下記の設定を行います。

Auto Analysis Workflow

自動解析・データ出力の設定 ラン後に自動的に解析データが出力されます。 出力するデータおよび保存先は任意に指定できます。

E-mail Notifications

ラン状況を登録してあるメールアドレスへ 自動送信する設定 7900HTシステムが施設のIT管理担当者によってネットワーク接続されている場合に

利用できます。

設定についてはネットワーク知識が必要なため、施設のIT管理担当者にご相談ください。

各設定を確認・変更する場合はBatch Settings ボタンをクリックします。 設定確認後、ダイアログ内のOK ボタンをクリックしてBatch Settings を閉じます。

<u>ランの開始</u>

 排出スタックにプレートが残っていないことを確認し、使用スタックにプレートをセットします。 セット方法はp.8をご参照ください。 画面右下のStart Batchボタンを押してランを開始します。

SDS Automation Control	ller			
the first internation that	190			
THE LAR ENVIRONMENT TODOL HEID				
	- W.			
Instrument Overview	Plate Queue Fun Status Processed Plates			
Satur: Ide	5 Rates Currently Queued			
Blad: Type: 304 Welk Blad:	# Barcode Plate Nome	Pun Type		
Serial Number: 200027	1 394N03850R Dr/Applied Borystener/SE6 Decumental/2.3Auto/394N00850P	R.ode Real Tree		
044.044	2 38493385PS Dr/yppled Booystens/(SE6 Documents/): 36453/38493085PS.	ada Real True		
baut scales	A Setting Compared Booydenergics Excurrency: autoralettal A Setting Has Disposed Booydenergics Excurrency: autoralettal	Loss Pearline		
Quinted Habits	6 3940038597 DrAppled Bosystens(505 Documents); 34403(3940036597	sis Peal Title		
Processed nates: 0				
Patrician Description				
Currently Running Plate				
Run Time Remoting:				
Name:				
Barcode:				
Run Type:				
	Taxilla Addition	I maxim I remaindered I mean at I		
	-Automation Robot Settings			
	Select the stacks that contain the plates for this batch			
	IZ Stack1 Stack2 Stack3 Stack4 Restack places after run			
	Chable batch settings			
	IZ Auto-Analysis Workflow IZ E-nail notifications	Eatch Settings		
		Open/Close Deck		
				1
			One of Class Deer	Chave Datab
			Open/Close Door	Start Batch
			Open/Close Door	Start Batch
			Open/Close Door	Start Batch
			Open/Close Door	Start Batch
			Open/Close Door	Start Batch
			Open/Close Door	Start Batch
			Open/Close Door	Start Batch
			Open/Close Door	Start Batch

2. ランを開始するとRun Status タブ内にラン中の温度モニター状況が表示され、登録されたドキュメントの 自動ランが行われます。途中で終了する場合は画面右下のStop Batch ボタンをクリックします。

e Edit Enstrument Tools Help		
: 🛯 🖓 🖷 🖉 🖻 🕷 💊	W	
strument Overview tabusi Plate Rum lock Type: 304 Wells Block end Manhare: 20027	Plate Queue Run Statur Processed Plates - Cycle Information Stage: 1 Repetitors: 1 Stage: 1 State: Hidding	indument langerature Sample: 50.0 Cover: 204.6
atch Stabue	Time: 01:41	Bod: 50.0
recessed Plates: 0		Contraction of the second s
Mall Microsoft Marcine (Marcine) (Ma		
	8.0	103 0.4 0.80 0.84 Time (th) : sm)
	Interval . E Standa . E Cover . E E Nova	

解析終了後のプレートドキュメント

ラン終了後のプレートドキュメントに蛍光データが書き込まれますので、個別に再解析を行うことが可能です。 なお、解析については納品説明補足資料・簡易操作ガイドもしくは各アッセイに対応した英文操作ガイドを ご参照ください。