

Applied Biosystems 7900HT Fast Real-Time PCR System

TaqMan[®] Low Density Array 簡易操作ガイド

SDS software version 2.3





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。

TaqMan Low Density Array:

This product is a Licensed Probe. Its use with an Authorized Core Kit and Authorized Thermal Cycler provides a license for the purchaser's own internal research under the 5' nuclease patents and basic PCR patents of Roche Molecular Systems, Inc. and F. Hoffmann-La Roche Ltd. No real-time apparatus or system patent rights or any other patent rights owned by Applera Corporation, and no rights for any other application, including any in vitro diagnostic application under patents owned by Roche Molecular Systems, Inc. and F. Hoffmann-La Roche Ltd claiming homogeneous or real-time amplification and detection methods, are conveyed expressly, by implication or by estoppel.

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.

TaqMan is a registered trademark of Roche Molecular Systems, Inc.

© 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またはパッケージインサート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

Relative Quantification ~LDAとRQソフトウェアによる検量線を用いない相対定量~

<u>概要</u>

TaqMan[®] Low Density Array (LDA)は2µ1の反応ボリュームでリアルタイムPCRを行うことができるシステムで、Comparative Ct 法を用いて遺伝子発現定量を行います。

LDAではランダムプライマーを用いての逆転写反応液を鋳型とした、2ステップRT-PCRで使用することが必要です。

また必ず各カード内に内在性コントロール遺伝子が設定されている必要があります。

1度に解析を行う組み合わせの中に、比較相対定量のためのキャリブレーターサンプルを準備します。RQソフトウェアでは1度に10枚のカードを解析することができます。

LDAでは次のステップで検討を行います。

cDNA合成

反応液の調整

サンプルの注入

カードの遠心

シーリング

SDS 2.3ソフトウェアのセットアップ/ラン開始

RQ Manager 1.2ソフトウェアによるデータ解析

<u> AACt (RQ) ドキュメントを用いたデータ作成の注意点</u>

LDAを用いた比較C_T法(ΔΔ C_T法)による相対定量実験を行う際、最初にΔΔ Ct(RQ)ドキュメントを 用いてリアルタイムPCRアッセイを行い、RQ Manager 1.2ソフトウェア上で解析を行う必要があります。

<u>比較 C_T 法($\Delta\Delta C_T$ 法)とは</u>

比較 C_T 法($\Delta\Delta C_T$ 法)とは、基準としたサンプルとの C_T 値(Threshold Cycle値)の差から 相対値を求める相対定量法です。

比較C_T法(△△ C_T法)を用いて相対定量を行うためには、下記の2つの条件を満たしている ことが必要です。

- PCR増幅効率がほぼ100%であること
- ターゲット遺伝子と内在性コントロール遺伝子の2つの反応系でPCR増幅効率が ほぼー致していること

*弊社推奨条件に従って設計されたプライマー・プローブセットは、上記条件を満たしている 可能性が非常に高いと考られます。

<u>比較C_T法(ΔΔ C_T法)と検量線法の違い</u>

比較C_T法(ΔΔ C_T法)と検量線法の違いは下表の通りになります。

	長所	短所
比較C _T 法(ΔΔ C _T 法)	検量線作成用サンプルの 必要がないので、未知サン プル処理数が増え、コスト が下がる。	PCR増幅効率が100%だと 仮定しており、実際のPCR 増幅効率が実験結果に反映 されない。
検量線法	毎回PCR増幅効率を反映 して検量線が作成される ため、得られた定量値の正 確性が高い。	毎回の実験毎に検量線を 作成することが必要なため、 未知サンプル処理数が減り、 コストが高くなる。

それぞれの長所と短所をご理解いただき、ご検討ください。

<u>比較C_T法(ΔΔ C_T法)に関する資料</u>

比較C_T法(△△ C_T法)の原理等の詳細については "ABI PRISM 7700 Sequence Detection Systems User Bulletin #2"をご参照ください。 この資料は弊社ホームページ(www.appliedbyosystems.co.jp)からダウンロードする ことができます。

当社へのお問い合わせ

アプリケーションサポートへのお問い合わせ お問い合わせ先の電話番号、電子メールアドレスは下記の通りです。 アプリケーションサポート フリーダイアル TEL: 0120-477-392 FAX: 0120-477-120 電子メール : jptechsupport@appliedbiosystems.com

<u>用語</u>

内在性コントロール(Endogenous Control)

相対発現定量では、テンプレート量に影響を与える核酸の抽出及びRT効率の補正のために、ターゲット遺 伝子の他に、内在性コントロール遺伝子の定量を行います。コントロール遺伝子は18s rRNA, β-actin, GAPDH などが用いられますが、対象の実験系で発現が変動しない遺伝子を用いることが必要です。

キャリブレーターサンプル

Comparative C_T 法による解析では、基準になるサンプルをキャリブレーターとして設定することで、そのサンプルに対する相対値として定量結果を算出します。タイムコースの実験では0時間サンプルもしくは刺激前サンプル、また組織特異性の実験では比較対照にしたい組織由来のサンプルなどをキャリブレーターサンプルに設定します。

レプリケート

統計学上精度の高い定量結果を得るために、複数のレプリケートを設定することを推奨します。

棄却検定による解析対象ウェルからの自動除去

SDSソフトウェアの機能で、レプリケート数を3以上回収した場合に、スミルノフ-グラブスの棄却検定アルゴリズムに基づき、シングルウェルではC_T値、マルチプレックスランでは△C_T値が外れている場合に、自動的に解析対象から外します。解析の設定で棄却検定を用いるかどうかを設定します。

<u>cDNA合成</u>

High Capacity cDNA Reverse Transcription Kitを用いてcDNA合成を合成します。 RNase Inhibitorなし P/N: 4368814・200反応, 4368813・1000反応 RNase Inhibitorあり P/N: 4374966・200反応, 4374967・1000反応 High Capacity cDNA Reverse Transcription Kit Protocol(P/N: 4375575)に準じて反応を 行ってください。

<u>反応液の調整</u>

1.1ポートにつき以下の容量で反応液を調整します。

コンポーネント	容量(uL)/リザーバー
cDNAサンプル*	a uL
RNase/DNase フリー水	(50-a) uL
TaqMan Universal Master Mix(2X)	50 uL
Total	100 uL

•RNA量換算で30~1000ngのcDNAを使用します(通常100~200ng)。

- 2. 1.をやさしく攪拌します。
- 3. スピンダウンします。

<u>サンプルの注入</u>

- 1. LDAのパッケージをあらかじめ室温に戻しておき、取り出します。
- 2. アルミ面を下にして実験台上に置きます。
- 3. 100 µ Lの反応調整液を Fill ポートからゆっくりと注入します。 (カードのフォーマットに従って、1検体あたり1~8ポートに注入します。)





<u>カードの遠心処理</u>

- 1. カードホルダーにLDAカードを挿入します。空いたポジションにはブランクカードを入れます。
- 2. "This side Out "が外側になるようにホルダーをバケットに入れます。



3. バケットを遠心機にバランスよくセットし、331g(室温)にて1分間遠心します。

Parameter	QUIKSet	EASYSet
Up ramp rate	3	9
Down ramp rate	N/A	9
Rotational speed	1200 rpm	1200 rpm
Centrifugation time	2 x 1 min (0.01 on display)	2 x 1 min (0.01 on display)

- 重要! 1200 rpmを超える回転数で遠心したり、3分以上遠心することは絶対に避けてください。 カードが変形、破損する恐れがあります。
- 4. 3が終わったら、気泡が入っていないか確認し、もう一度331g(室温)にて1分間遠心します。

シーリング

- 1. シーラーを水平の台上に置きます。
- 2. スタートポジションが手前(自身に近い側)に来るように置きます。
- 3. Carriage を Starting ポジションにセットします。



- 4. カードをホイルの部分が上になるようにシーラーに置きます。
- 5. Fillポートが向こう側になるようにシーラー上に置き、やさしく押してPinでカードが固定されるようにします。



6. Carriageを向こう側にゆっくり押します。

(土台に手をかけるとCarriageにはさまれることがありますので注意してください)



- カードに8本の窪みができ、シールされます。
 (カードがセットされた状態でCarriageをEnd ポジションから動かさないでください)
- 8. カードを取り出します。
- 9. Fillリザーバーの部分をはさみで切り取ります。



<u>新規プレートドキュメントの作成</u>

1. デスクトップ上のショートカットをダブルクリックし、SDS Software を起動します。



- 2. File メニューより New を選択します。
- 3. New Document が表示されます。

Assay:		-(4
ontainer:	384 Wells Taqman Low Density Array	-(
emplate:	Blank Template	
	Browse	
arcode:		

- 4. Assay のドロップダウンリストより ΔΔCt (RQ) を選択します。
- Container のドロップダウンリストより、反応を行うプレートタイプを選択します。
 384 wells TaqMan Low Density Array を選択します。
- 6. **OK** をクリックします。
- 7. 新規のプレートドキュメントが開きます。

プレートドキュメントについて

プレートドキュメントは大きく3つのフィールドに分かれています。



- A) グリッド:ウェルを選択するフィールドです。
 このフィールドで選択したウェルに対して設定を行うことができます。
 また、このフィールドで選択をしたウェルのデータを表示させることができます。
- B) システムテーブル:ウェル毎の設定や解析結果を確認するフィールドです。 各ウェル毎の設定や、ラン終了後には定量値等がここに表示されます。
- C) タブを切り替えて作業を行うフィールドです。
 Setupタブ: Detectorの登録と、各ウェルに対する設定を行います。
 Instrumentタブ: サーマルサイクラー条件の設定を行います。

 SDS Setupファイルをインポートします。 Setup ファイルはLDA上にアプライされている遺伝子検出用プライマー/プローブセット(Detector)の 位置情報が記録されたファイルになります。 このファイルを用いることで、プレートドキュメントの設定を非常に簡単に行うことができます。 ファイル名は [SDS_Product Number.txt] という形式になっています。Product Numberはご注文ごとに 割り振られる固有の番号です。



- a) LDAに付属しているCDをセットし、SDS setupファイルを確認します。
- b) Fileメニューより Import を選びます。Importダイアログボックスが表示されます。
- c) CDを開いてSetupファイルを選択し、Import をクリックします。
- Well gridより全てのウェルを選択します。
 必要に応じてTaskをTarget 又は Endogenous Control に変更してください。

Setup	Instrument			
Wel(s)	: A1-P24			?
Sample	Name: * Mixed *			_
Use	Debactor	Reporter	Task	Color
F	4342376-GAPDH	FAM	Endogenous Control	
F	CD44-Hs00174139_m1	FAM	Target	
	CD63-Hs00168+66_m1	FAM	Target	× .
F	CD86-Hs00199349_m1	FAM	Tarpet	
F	COL1A2-Hs00164099_m1	FAM	Endogenous Control	
F	COL6A1-Hs00242448_m1	FAM	larget	
F	CSF2-Hs00171266_m1	FAM	Target	
F	EGR1-Hs00152928_m1	FAM	Target	
F	FGF2-Hs00256645 m1	FAM	Target	

10. Well grid で同一サンプルが入っているウェルを選択し、サンプル名を入力します。

サーマルサイクリング条件の設定とランの開始

1. Instrument タブをクリックすると LDA専用の温度条件が表示されます。

AmpErase UNG	ApmliTaq Gold Each of		40 Cycles
Activation	Polymerase Activation	Melt	Anneal/Extend
50°C ∕ 2min 100% ramp	94.5℃/10min 100% ramp	97°C∕30sec 50% ramp	59.7°C ∕ 1min 100% ramp

※温度、時間、反応容量等、サーマルサイクリング条件はこのままで使用します。

- ランを行う前にドキュメントを保存します。
 ランをするために必要なプレートドキュメント保存形式は拡張子.sdsとなります。
 テンプレートファイルとして保存する場合は拡張子.sdtを選択して保存します。
 - 重要 テンプレートファイルは必ずランを行う前のプレートドキュメントから作成して下さい。 ラン後の.sdsファイルからテンプレートファイル(.sdt)を作成しますと不完全なファイルとなり、 ランが正常に行われませんのでご注意ください。
- Instrumentタブの Open/Close ボタンをクリックし、トレイを移動させます。
 A1ポジションが左上、ノッチコーナーが右上、バーコードが手前から見えるようにカードの アルミ面を下にしてトレイ上に載せます。



- 4. Start ボタンをクリックして、ランを開始します。
- 5. ランが終了したら、Save をします。

RQ Manager 1.2 ソフトウェアを用いたRelative Quantification Study

概要

△Δ Ct (RQ)ドキュメントをRQ managerソフトウェア上で解析することで、比較C_T法(ΔΔC_T法)を 用いた相対定量値の算出を行うことができます。

RQ Manager ソフトウェアでは最大10枚の△△ Ct (RQ)ドキュメントを読み込み、同時に解析を 行うことができます。

重要

複数枚の △△ Ct (RQ)ドキュメントを同時に解析する場合、全てのドキュメントで同じサーマルサイクリン グ条件(ステップ、サイクル数、反応容量、9600 Emulation の有無)で設定することが必要です。 異なる設定があるとRQ Studyドキュメントに同時に読み込んで解析を行うことができませんのでご注意く ださい。

RQ Study ドキュメントの新規作成

1. デスクトップ上のショートカットをダブルクリックし、RQ Manager Software を起動します。



- 2. File メニューより New Study /From SDS File を選択します。
- 3. New Study ダイアログが表示されます。
 - a) Container Type のドロップダウンより、384 Wells TaqMan Low Density Array を選択します。
 - b) この RQ Study のファイル名を入力します。
 - c) Add Plate(s) をクリックします。

Container Type	384 Wells Taqm	an Low Density Array		
Study Name	Untitled			- (3b)
ates:				
Nar	ne	Runtime	Туре	T
ⅆ Plate(s)		- 30		
Add Plate(s).]	30		

- 4 Add Plates ダイアログボックスが表示されます。
 - a) 画面左側にコンピュータ内に保存されているファイルが表示されます。解析する∆∆ Ct(RQ) ドキュメントを選択します。
- b) Add ボタンをクリックします。
- c) OK ボタンをクリックし、ダイアログボックスを閉じます。

File Search	Chilleniad Riscustere/CDC Decumenta/Comple RO
S My Computer	Add >> C. Applied biosystems (SUS Documents (Sample RG
🗄 🔩 31/2 Floppy (A:)	
E Secol Disk (C:)	Remove
🗄 🧰 ABI	
Applied Biosystems	4b
🖶 🧰 SDS Documents	
Appliedbiosystems	
Decc Deccc Deccc Decc Deccc Decc Decc Decc Decc	
B Columnation of the octuality of the second s	Evicting Distan In The Study
Program Files	Existing Plates in the Study
B C RECYCLER	
B C rho inst	
I C Documents	
B C SDS 1	
III 🛅 si659	
System Volume Information	
I Contemp	
I CONTRACTOR INTERNAL	
🗄 🥝 DVD Drive (D:)	
🗷 🎱 CD-RW Drive (E:)	
My Network Places	
Documents	(4c)
B Share-to-Web Upload Fobler	
4352716B_mtr	
Number of Distan Obuly and holds 40, 0 feet	
Number of Prates Study can hold: 10 - 9 free	OK
—	

- 5. New Study ダイアログボックスにドキュメントが取り込まれていることを確認し、OK をクリックします。
 - 新規RQ Study ドキュメントが開きます。

取り込める残り枚数が表示されます。

RQ Study ドキュメントでの解析

1. Analysis メニューから Analyze Settings を選択するか **国**をクリックして解析パラメータの設定を行います。 Analysis settingには3つのタブ画面があります。

Studyタブ

Γ Omit Replicates whose Δ0	1 <= 1.0 (multi	olex only)			
Gene Expression Settings for Stu	idy				1a
Endogenous Control Detector:	GAPDH		-		-
Cont	rol Type: C Multiple:	ed 🕼 Non-r	nultiplexed	1	
RQ Min/Max Confidence:	99.0%		7	—(1
					-
Flag Condition and Omit Settings					12002
Flag Condition and Omit Settings Flag Condition and On	nt Studies When	Flag	Condition	Flag Conditi	Omit
Flag Condition and Omit Settings Flag Condition and On Nell is an outlier (GBO)	nit Studies When	Flag	Condition	Flag Conditi	Omit
Flag Condition and Omit Settings Flag Condition and On Well is an outlier (GBO)	nit Studies When	Flag	Condition	Flag Conditi	0

- a) 内在性コントロールの指定を行います。ドロップダウンリストの中から任意に選択できます。
- b) データばらつきの統計学的な信頼度を選択します。
- c) フラッグや自動データ除外設定を行います。

Detector タブ

Ct Analysis	-				
Manual Ct	Threshold:				
Automatic Basel	ine				
C Manual Baseline	Start (Cycle):	3	End (Cycle): 15 ÷	

d) ベースラインおよびスレッショルドラインの設定を行うタブになります。Automatic Ct を 選択すると、全 Detector のベースラインとスレッショルドラインが自動決定されます。

Plate タブ

自動除外やQCチェックのパラメータ設定タブになります。 具体的なフラッグの内容については英文ガイド["]Maintenance and Troubleshooting Guide["] p.126からをご参照ください。

2. すべての設定を行ったのちOK もしくは Apply をクリックすることで解析が行われます。

 Analysis メニューから Analysis All を選択するか ・ をクリックして解析を行います。

 (2.のステップですでに解析が行われている場合はボタンはグレイアウトします。)
 解析が終了すると、下図のような画面が表示されます。



A) RQ Detector Grid

ここにはRQ Study ドキュメントに設定された Detector が表示され、クリックするとその結果が C)の画面に表示されます。

B) RQ Sample Grid

A) で選択した Detector で設定されているサンプル及び △△ Ct 法による定量結果その他の情報が 表示されます。

C) RQ Results Panel

A) で選択した Detector の Amplification Plot や Gene Expression のグラフが表示されます。

 A) のグリッドで Detector を一つずつ選択し、C) に表示される増幅曲線およびスレッショルドラインが 適切な位置に設定されているか確認して下さい。もし適切でない位置に設定されている Detector が 存在していた場合、p.15の Analysis Settings 画面にて Manual Ct に切り替えて再度解析してください。 ベースラインおよびスレッショルドラインの設定についてはp.17, 18をご参照ください。







[短すぎるベースライン] ベースラインのEnd cycle を大幅に 越えてから増幅曲線が始まって いるので、End Cycle 値を 増やす必要があります。



[長すぎるベースライン] 増幅曲線がベースラインのEnd cycle以前に始まっているので、 End Cycle 値を減らす必要が あります。

―スレッショルドラインの設定―

[正しく設定されたスレッショルド] 設定されたスレッショルドは、 増幅曲線の指数関数増幅領域内に あります。 スレッショルドが最適値を上回った り下回って設定されると、複製グル ープの標準偏差が増加します。



[低すぎるスレッショルド]

設定されたスレッショルドは、増幅曲線の 指数関数増幅領域より低い領域にあります。 この場合、スレッショルドが正しく設定 された場合のプロットよりも、はるかに 多くの標準偏差が発生します。 スレッショルドバーを指数関数増幅 領域内に引き上げてください。



[高すぎるスレッショルド]

設定されたスレッショルドは、増幅曲線 の幾何学的相より高い領域にあります。 この場合、スレッショルドが正しく設定 された場合のプロットよりも、はるか に多くの標準偏差が発生します。 スレッショルドバーを指数関数増幅 領域内に下げてください。



- 5. Gene Expression タブでの相対発現定量結果の確認 △ △ Ct 法による相対発現定量結果が表示されます。縦軸にキャリブレーターサンプルに対する 相対量のlog 値、横軸に Detector やサンプルが表示されます。
 - a) Gene Expression タブに切り替えます。
 - b) 表示させたい Detector を選択します。
 - c) グラフの表示形式や Calibrator サンプルを変更する場合にドロップダウンリストから選択します。



6. グラフのスケールは、グラフ軸上でダブルクリックするか Analysis メニュー内の Graph Settings を 選択することで表示されるダイアログボックス上で変更可能です。

aph Settings			
-X Axis -X Axis Auto So Bar Width:	ale 12 pixel	s	
Y Axis	I⊽ Auto Sca Minimum	le -5,0	
	Maximum	5.0	
Save Settings /	As My Default	ок	Cancel

相対定量の測定結果の表示は以下の意味を持ちます。

①「相対発現量バー」+「エラーバー」による表示



キャリブレーターサンプル及びテストサンプルのターゲット遺伝子及びコントロール遺 伝子において全ての有効なC_T値が算出されている(設定したC_T値の最大値より小さい) 場合、相対発現量+エラーバー(信頼水準より算出した相対量のとりうる最大/最小値の 範囲)で表示されます。

②「相対発現量の最大値バー」+「↓」による表示



キャリブレーターサンプルのターゲット遺伝子及びコントロール遺伝子は有効なC_T値 が算出され、かつテストサンプルのターゲット遺伝子のC_Tが算出されていない場合、とり うる「最大レベルの相対発現量」+「↓」で表示されます。このことは、測定を行うテストサ ンプル中に正確な相対定量を行うために必要なターゲットcDNAが含まれていないことを 示しており、テストサンプルのテンプレート量を増やして再実験する等が必要になります。

③「相対発現量の最小値バー」+「↑」による表示



テストサンプルのターゲット遺伝子及びコントロール遺伝子は有効なC_T値が算出され、 かつキャリブレーターサンプルのターゲット遺伝子のC_Tが算出されていない場合、とりう る「最小レベルの相対発現量」+「↑」で表示されます。このことは、キャリブレーターサン プル中に正確な相対定量を行うために必要なターゲットcDNAが含まれていないことを示 しており、キャリブレーターサンプルを変更して再解析を行うか、キャリブレーターサンプ ルのテンプレート量を増やして再実験する等が必要になります。

その他の結果の場合には、相対量の表示は行われません。cDNA量の変更や実験系の検討が必要になります。

<u>解析結果の出力</u>

- 1. File メニューの Export から Results を選択して、下記の中から出力する データを選択します。
 - Amplification data
 - Expression data
 - Clip data
 - •Results data
- 2. 出力ファイルのファイル名を入力し、任意の場所を選択し、Save ボタンを クリックします。
- 3. タブ区切りのテキストファイルとして出力されます。

アプライドバイオシステムズジャパン株式会社