

# SureCycler 8800

# 日本語クイックガイド

バージョン A, 2011 年 5 月



Agilent SureCycler 8800 ユーザーガイド

# **Agilent Technologies SureCycler 8800 Installation Instructions**

1 広いスペースで SureCycler の外箱を開けてくだ さい。 一番上に動作テスト合格証明書 (Functional Test Certificate)、CDおよびイン ストールポスターが入ったアクセサリートレイ があるのでそれを取り出します。 2 本体の両脇から手を入れ、必ず本体に手をかけて 持ち上げてください。梱包材ごと本体を持ち上 げ、箱から出します。



4 サーマルブロックの箱を開けます。(A) サーマルブロック上部にある梱包材を取り除きます。
 (B) ヒートシンクを折らないように気をつけて、サーマルブロックを箱から取り出します。





a form to be a second	and the second	and property of
inte Nutera	201000000000000000000000000000000000000	

💥 Agilent Technologies











The Measure of Confidence

Agilent Technologies

www.agilent.com/genomics/8800

Agilent SureCycler 8800 ユーザーガイド

# 注意

© Agilent Technologies, Inc. 2011 本ユーザーガイドは米国著作権法 および国際著作権法によって保護 されており、Agilent Technologies, Inc. の書面による事前の許可なく、 本書の一部または全部を複製する ことはいかなる形式や方法(電子 媒体による保存や読み出し、外国 語への翻訳なども含む)において も、禁止されています。

#### テクニカルサポート

テクニカルサポートについては Agilent

<u>email japan@agilent.com</u> までご 連絡ください。連絡先情報は14ペ ージに記載されています。

#### 保証

このマニュアルに含まれる内容は、 「現状のまま」提供されるもので、 将来の改定において予告なく変更 されることがあります。また、 Agilent は、適用される法律によっ て最大限に許可される範囲におい て、このマニュアルおよびそれに 含まれる情報に関して、商品性お よび特定の目的に対する適合性の 暗黙の保証を含み、それに限定さ れないすべての保証を明示的か暗 黙的かを問わず一切いたしません。 Agilent は、このマニュアルまたは それに含まれる情報の所有、使用、 または実行に付随する過誤、また は偶然的または間接的な損害に対 する責任を一切負わないものとし ます。Agilent とお客様の間に書面 による別の契約があり、このマニ ュアルの内容に対する保証条項が この文書の条項と矛盾する場合は、 別の契約の保証条項が適用されま す。

#### 技術ライセンス

このマニュアルで説明されている ハードウェアおよびソフトウェア はライセンスに基づいて提供され、 そのライセンスの条項にしたがっ て使用またはコピーできます。

#### 安全に関する注意

#### 注意

注意は、危険を表します。これは、 正しく実行しなかったり、指示を 順守しないと、製品の損害または 重要なデータの損失にいたるおそ れがある操作手順や行為に対する 注意を喚起します。指示された条 件を十分に理解し、条件が満たさ れるまで、注意を無視して先に進 んではなりません。

#### 警告

警告は、危険を表します。これは、 正しく実行しなかったり、指示を 順守しないと、人身への障害また は死亡にいたるおそれがある操作 手順や行為に対する注意を喚起し ます。指示された条件を十分に理 解し、条件が満たされるまで、警 告を無視して先に進んではなりま せん。

# このユーザーガイドでは…

本ユーザーガイドはAgilent SureCycler 8800の設置方法および使用方法について説明しています。

- 1、始める前に この章では、使用を開始する前に読む必要のある情報を記載しています。
- 2、インストール

この章では、SureCycler 8800の設置方法について説明しています。

3、基本操作

この章では、サンプルの準備・ロードの方法、ソフトウェアの操作方法、ユーザーアカウント の設定などについて説明しています。

4、PCR プログラム

この章では、PCR プログラムの作成、実行、インキュベーションの実行、レポートの表示など について説明しています。トラブルシューティングもこの章に含まれています。

5、メンテナンスとソフトウェアの更新 この章では、機器のメンテナンスおよびソフトウェア更新のアップロードに関する方法につい て説明しています。

SureCycler 8800 日本語クイックガイド 3

# 目次

1. <b>始める前に</b>	9
梱包の内容	10
安全に関する注意事項	10
インジケーターLED ライト 13	
概要	13
SureCycler8800 技術サポート	14

# 2. インストール

インストール		15
ステップ 1	機器の設置場所の選択	16
ステップ 2	機器の開梱	16
ステップ3	サーマルブロックユニットの本体への取り付け	16
ステップ4	電源への接続	17
ステップ 5	オプションデバイスの接続	17
ステップ6	電源投入	17
ステップ7	ネットワーク接続(オプション)	17
サーマルブロ	ックの交換	18
サーマルブ	ロックの外し方	18

リーマルノロックの外しり	
サーマルブロックの取り付け方法	18

3.	基本操作	19
	サンプルの準備とロード	20
	サンプルの準備	20
	サンプルのロード	21
	Agilent プラスチック製品および消耗品	21

ソフトウェア操作	22
タッチスクリーン操作	22
ホーム画面	22
ユーザーアカウント	24
アクセスレベル	24
ユーザーアカウントへのログイン	25
ユーザーアカウントの管理	26
ネットワーク設定とリモートアクセス	29
装置の IP アドレスの確認・編集	29
別の装置からのアクセス	30
PC から装置へのアクセス	31
装置本体の設定	32
ホットトップの設定	32
サーマルサイクラーの名前	33
日付と時刻	33

### 4. PCR プログラム

新規 PCR プログラムの作成	36
手動でプログラムを作成するには	36
ウィザードを使用して PCR プログラムを作成するには	39
384 ウェル用に PCR プログラムを作成する際のヒント	41

コマンドの説明	42
Hot Start (ホットスタート)	42
Cycle Wizard(サイクルウィザード)	43
Start Cycle(サイクル開始)	44
Temperature Step (温度ステップ)	45
End Cycle(サイクルの終了)	46
Gradient Step (グラジエントステップ)	45
Touchdown(タッチダウン)	48
Storage/Pause Step(保存・一時停止ステップ)	50

プリインストールされたプログラムの使用	51
プリインストールされたプログラムの利用	52
プリインストールされたプログラムの編集	52
PCR プログラムの編集	53
プログラムのパラメータを編集するには	53
PCR ステップをコピーするには	54
PCR ステップを追加または削除するには	54
PCR プログラムを削除するには	54
PCR プログラムの実行	55
PCR プログラムを実行するには	55
実行中のプログラムを中止または一時停止するには	55
インキュベーションの実行	56
インキュベーションの設定と実行	56
レポートの表示	57
レポートを表示するには	57
レポートを削除するには	57
トラブルシューティング	58
スタンダード・ランプレートを使って、	
少量で反応させた際に PCR 産物の収量が低い場合	58
プログラムの実行中にブロック温度が表示されない場合	58
プログラムの実行が始まらない場合	58
チューブの内部に結露があった場合	59
シールしたマイクロプレートでサンプル液量が減少した場合	59
メモリスティック、キーボードなどの USB デバイスが認識されない場合	59
装置の電源が入らない場合	59
画面が点灯せず、何も表示されない場合	60
画面が点灯するものの、何も表示されない場合	60

5. メンテナンスとソフトウェアの更新	61
メンテナンス	62
クリーニング	62
ヒューズの交換	62
ソフトウェアの更新	63

SureCycler 8800 日本語クイックガイド 8



1. 始める前に

梱包の内容	10
安全に関する注意事項	10
インジケーターLED ライト	13
概要	13
SureCycler 8800 技術サポート	14

この章では、本機の使用を開始する前に読む必要のある情報を記載しています。 詳細については、16ページの「インストール」を参照してください。

# 梱包の内容

表 1 梱包の内容

部品	数量
SureCycler 8800 機器ベース	1
サーマルブロックユニット(96 ウェルまたは 384 ウェル )*	1
電源コード	1
機能テスト証明書	1
クイックセットアップポスター	1
設定およびユーザーガイド CD	1

\*サーマルブロックユニットは部品番号 G8810A(96 ウェルブロック)および G8820A(384 ウェルブロック)で 追加購入可能

# 安全に関する注意事項

電気

以下を含めた標準的な電気的安全に関する注意事項を守る必要があります。

・必要に応じてすぐに電源を外すことができるような場所に機器を設置するようにしてください。

- ・最初に機器に電源を入れる時には、電圧が適切なレベルであることを確認する必要があります。
- ・デバイスは接地された電源コンセントに接続する必要があります。
- ・濡れた手でスイッチやコンセントに触れないでください。
- ・電源コードを外す前に機器をオフにしてください。

・こぼれた液体等を清掃する場合や、内部コンポーネントを修理したり、サーマルブロックを交換する場合には、 機器の電源を抜いてください。

- ・接地されていない電源に接続しての使用はしないでください。
- ・電気コンポーネントを修理するには資格が必要です。

#### 1. 始める前に

#### 安全に関する注意事項

#### 液体および試薬

・サイクラーの外側にあるサーマルブロックに液体を充填する際には液体が機器に浸透しないようにしてください。

・爆発性、可燃性、および反応性の物質を機器でインキュベーションまたはサーマルサイクルにかけないでくだ さい。

・病原性物質や放射性物質、その他健康に有害な物質を処理するときには関連する安全規制を順守する必要があります。

・機器を液体に浸さないでください。

#### やけどの危険

・サーマルブロックやホットトップには触れないでください。これらのエリアは急速に 50°C 以上の温度になります。温度が 30°C 以下になるまで、ホットトップを開かないでください。

・十分な温度安定性がない材質の製品 (プレート、プレートシール、フォイル、マット)を使用しないでください。(最大で 120°C 程度になります)

#### 動作環境

・デバイスの通気スロットは常時通期可能な状態に保つ必要があります。機器の周囲に少なくとも 10 cm のス ペースを空けてください。

・機器の周囲温度は 20 ~ 30°C、湿度レベルは 20 ~ 80 %で結露しないようにして下さい。

・有毒または爆発の危険性がある環境で機器を操作しないでください。

#### 機器定格

- 汚染度 2
- ・設置カテゴリ 2
- ・高度 2000 m
- ・湿度 20 ~ 80%、ただし結露しないこと
- ・電源 100 ~ 240 V、50/60 Hz、11 A
- ・温度 20 ~ 30°C
- ・室内使用専用

#### 1. 始める前に

#### 安全に関する注意事項

#### 静電気放電

SureCycler 8800 は静電気の影響を受けやすい機器です。8000 ボルト以上の静電気放電は機器の USB ポートの 正常な動作に影響を及ぼす恐れがあります。高い静電気放電があるような環境で作業をする場合には注意が必要 です。高い静電気放電がある環境では、接地されたリスト ストラップの着用や、他の静電気防止策を取ったう えで機器に触れる必要があります。ESD STM5.1-1998 クラス 3B.

#### 安全記号

以下の電気 / 安全記号は機器に表示されているものです。



# インジケーター LED ライト

機器の前面にはステータスインジケーター LED があります (右上隅)。この LED ライトのステータスコードを 以下の表に示します。

表 2 ステータスインジケーター LED の表示

表示	機器ステータス
オフ	サイクラーはアイドル状態です。
緑色に点滅	サイクラーはプログラムまたはインキュベーションを実行しています。
緑色点灯	サイクラーは一時停止しています。
	サイクラーがエラーを検出しました。ディスプレイでエラーメッセージをチェックし
	さらに詳細を確認してください。

### 概要

SureCycler 8800 はサーマルサイクラーであり、DNA 鋳型の増幅を行うポリメラーゼ連鎖反応 (PCR) および関 連したメソッドを実行するように設計されています。SureCycler 8800 は、サイクルごとの時間・温度の増減 (Cycle increment)、タッチダウン PCR、および温度勾配を含めた、きわめて複雑なサイクルプログラムを実行 することができます。カラー・タッチスクリーンで操作を行うソフトウェアインターフェースはインテリジェン トな Program Wizard を備えています。Program Wizard はプライマーおよび鋳型の情報から PCR プロトコルを 自動的に作成することができます。

#### 1. 始める前に

SureCycler 8800 技術サポート

# SureCycler 8800 技術サポート

電子メール : email\_japan@agilent.com

ホームページ www.genomics.agilent.com

電話:0120-477-111



2. インストール

インストール		15
ステップ 1	機器の設置場所の選択	16
ステップ 2	機器の開梱	16
ステップ3	サーマルブロックユニットの本体への取り付け	16
ステップ4	電源への接続	17
ステップ5	オプションデバイスの接続	17
ステップ6	電源投入	17
ステップ7	ネットワーク接続(オプション)	17
サーマルブロックの交換		

サーマルブロックの外し方	18
サーマルブロックの取り付け方法	18

この章では、SureCycler 8800の設置方法について説明しています

インストール
 インストール

# インストール

### ステップ1 機器の設置場所の選択

本機は水平できれいな場所に設置してください。以下の点にご留意ください。

- ・本機が安定した場所に設置されていること
- ・背面部には適当な間隔があり、通気が保たれていること
- ・本機は必ず壁または隣接する機器から少なくとも10cm以上離すこと
- ・本機は温度が20~30℃、湿度が20~80%で結露がない場所に設置されていること
- 本機周辺に爆発性あるいは可燃性のものがないこと

### ステップ2 機器の開梱

輸送用梱包箱を開き、本機および付属品を取り出します。 詳しくは、添付の設置ガイドをご参照ください。

### ステップ3 サーマルブロックユニットの本体への取り付け

1. サーマルブロックユニットの両側を両手でつかみ、装置本体の上に乗せます。

2. サーマルブロックユニットの両側に上から圧力をかけ、カチッとロックされるまで押し込みます。左右それ ぞれが確実にロックされるように、両側を押し込むようにしてください。

サーマルブロックユニットは、本体に正しい向きでセットされた場合にのみ所定の位置にロックされます。 サーマルブロックユニットを交換する際は、18 ページの「サーマルブロックユニットを交換する」を参照してく ださい。

# ステップ4 電源への接続

本機は接地付きコンセントに接続してください。

- 1. 装置本体の電源スイッチがオフになっていることを確認します。
- 2. 電源コードを機器背面にある電源コネクタに接続します。
- 3. 電源コードをコンセントに差し込みます。

# ステップ5 オプションデバイスの接続

必要に応じて、装置本体の前面にある USB ポートを使用してマウス、キーボード、USB メモリスティックなどを 接続することができます。

### ステップ6 電源投入

1.装置本体の左下にある電源ボタンを押し込みます

本機はいつでも電源を切ることができます。電源を切る前にソフトウェアを終了するなどの必要はありません。

### ステップ7 ネットワーク接続(オプション)

1. イーサーネットケーブルを装置本体の背面にあるイーサーネットポートに差し込みます。

2. イーサーネットケーブルの他方の端をネットワークポートに差し込みます。

# サーマルブロックの交換

SureCycler 8800 では、96 ウェルブロックと 384 ウェルブロックを交換して使用することができます。

### サーマルブロックの外し方

1. 装置本体前面の電源スイッチで電源を切ったのちに、電源ケーブルをコンセントから抜きます。

2. 本機上部にある灰色のボタンを押し、ふたを開けます。

3. 両手で左右からサーマルブロックユニットをつかみ、左上隅、右上隅にあるボタンをそれぞれの親指で押し込みます。

4. ボタンを強く押しながらサーマルブロックユニットを持ち上げ、装置本体から外します。

サーマルブロックユニットを取り外す際に、ふたが手前に倒れて手が挟まれないようご注意ください。サーマルブロックユニットを少し奥に倒しながら持ち上げることで、ふたが手前に倒れにくく なります。

### サーマルブロックの取り付け方法

1. サーマルブロックユニットの両側を両手でつかみ、装置本体の上に乗せます。

2. サーマルブロックユニットの両側に上から圧力をかけ、カチッとロックされるまで押し込みま

す。左右それぞれが確実にロックされるように、両側を押し込むようにしてください。

サーマルブロックユニットは、本体に正しい向きでセットされた場合にのみ所定の位置にロックされます。

3. 取り付けが完了したら再び電源に接続します。



3. 基本操作

サンプルの準備とロード	20
サンプルの準備	20
サンプルのロード	21
Agilent プラスチック製品および消耗品	21
ソフトウェア操作	22
タッチスクリーン操作	22
ホーム画面	22
ユーザーアカウント	24
アクセスレベル	24
ユーザーアカウントへのログイン	25
ユーザーアカウントの管理	26
ネットワーク設定とリモートアクセス	29
装置の IP アドレスの確認・編集	29
別の装置からのアクセス	30
PC から装置へのアクセス	31
装置本体の設定	32
ホットトップの設定	32
サーマルサイクラーの名前	33
日付と時刻	33

この章では、サンプルの準備・ロードの方法、ソフトウェアの操作方法、ユーザーアカウントの設定などに ついて説明しています。 サンプルの準備とロード

# サンプルの準備とロード

サーマルブロックおよびホットトップの損傷を防ぐために、耐熱性のあるチューブ、プレートをご使用 ください。不適切なチューブ、プレートを使用すると運転中に装置が損傷する可能性があります。推奨 のプラスチック製品・消耗品については 21 ページの「Agilent プラスチック製品および消耗品」をご参 照ください。

警告

やけどの危険:サーマルブロック、ホットトップ、チューブ、プレートは100℃近い温度に

達する場合があります。30℃以下になるまで手を触れないようにしてください。

### サンプルの準備

サンプルを準備する際には、以下のガイドラインに従ってください。

・プレートまたはチューブをサーマルブロックにロードする前に、プレートまたはチューブのキャップ
 を閉めてください。

・良好な結果を得るためには、サンプルをサーマルブロックにロードする直前に遠心機で短時間スピン ダウンしてください。

・プレートとシーリングフィルムの組み合わせで使用する場合には、反応中にコンプレッション・マットをシーリングフィルム上に載せてホットトップとの間に挟むことで、シーリングフィルムがプレート に圧着するようにしてください。コンプレッション・マットの注文に関しては 21 ページの「Agilent プラスチック製品および消耗品」をご参照ください。 本機はブロックタイプに応じて 96 ウェルプレート、384 ウェルプレート、シングルタイプの PCR チュー ブ、ストリップチューブが使用可能です。

- 1. 機器の上部にあるボタンを強く押し、サーマルブロックのふたを持ち上げます。
- 2. プレートまたはチューブをサーマルブロックにセットします。
- 3. ふたを閉め、カチッと音がするまで押し込みます。

# Agilent プラスチック製品および消耗品

以下の PCR プレート、チューブ、シーリングフィルムは SureCycler 8800 との互換性があります。

Agilent 製品番号	説明
401333	96 ウェルプレート 25 枚入り
410188	384 ウェルプレート 50 枚入り
410186	粘着シーリングフィルム、100 枚入り
410187	コンプレッション・マット、10 枚入り
410082	ストリップチューブ、200 uL,12 連チューブ 80 個入り
410086	ストリップキャップ、12 連キャップ(80 個入り

ソフトウェアの操作

# ソフトウェア操作

### タッチスクリーン操作

本機はカラー・タッチスクリーンが搭載されており、画面に表示されたボタンをタッチすることでソフ トウェアを操作することができます。USB 端子でマウスが接続されている場合には、クリックで操作す ることができます。以下の説明では、「タッチスクリーンを指でタッチする」および「USB 接続されたマ ウスでクリックする」操作を「タッチ」と表記します。プライマー配列、温度などのデータを入力する 必要がある場合には、タッチスクリーン上に仮想キーボードが表示されます。また、USB 接続のキーボ ードからデータを入力することも可能です。

### ホーム画面

本機の電源を入れるとタッチスクリーン上にはホーム画面が表示されます。このホーム画面は、装置本 体の制御やファイル操作、アカウント管理の起点となります。 以下にホーム画面の各ナビゲーションボタンの機能を記載します。

表3 ホーム画面ボタンの説明

ボタン	説明
New Program	新規 PCR プログラムを作成します。
Program	ファイルディレクトリを開き、保存済みの PCR プログラムを参照します。
	プログラムの編集、実行が可能です。
PCR Wizard	ウィザード機能を使用して新規 PCR プログラムを作成します。
Network Control	複数の機器にネットワークを介して接続します。
Incubate	ー定温度でのインキュベーションを行います。
Reports	レポートディレクトリを開き、保存されたレポートファイルを参照
	します。
Login	ログイン画面を開きます。
Settings	設定の変更、ソフトウェアのバージョンアップ、ユーザーアカウント
	管理を行います。

23 SureCycler 8800 日本語クイックガイド

ユーザーアカウント

### アクセスレベル

本機の電源を入れると、自動的に「Guest user」アカウントでログインした状態になります。

ユーザーアカウントとして、Administrator (管理者)、Registered user (登録ユーザー)および Guest user (ゲストユーザー)の3種類があり、それぞれ異なるレベルのユーザー権限があります。新規ユー ザーの登録、ユーザー権限やパスワードの設定・変更に関しては Administrator (管理者)のみが行う ことができます。

表 4	ア	ク	セス	レベ	ル
		-			

アクセスレベル	使用できる機能
Guest	・ゲストフォルダへのアクセス権限
	・ゲストフォルダに保存されているプログラムの作成、コピー、編集、
	および実行
User	・ゲストおよび個人フォルダへのアクセス権限
	・ゲストおよび個人フォルダに保存されているプログラムの作成、
	コピー、編集、および実行
Administrator	・すべてのゲストおよびユーザーフォルダへのアクセス権限
	・ユーザーアカウントのセットアップとメンテナンス
	・装置のセットアップとメンテナンス
	・すべてのレポートのメンテナンス、アーカイブ化

ユーザーアカウントへのログイン

#### Guest としてのログイン

Guest としてログインする場合にはパスワードは不要です。デフォルトでは、電源を入れるとゲストア カウントでログインする設定になっています。

Guest としてログインするとユーザー権限が制限され、誰でもアクセス可能なゲストフォルダを使用す ることになります。登録ユーザーとして装置を使用するためには、あらかじめ Administrator が新規ユ ーザーアカウントを設定する必要があります。登録ユーザーは、パスワードで保護された個人フォルダ を使用して PCR プログラムの管理をすることができます。

#### Administrator としてログインするには

1. ホーム画面で[Login (ログイン)]をタッチします。

Username ドロップダウンリストで Administrator を選択し、パスワードを入力します。デフォルトの Administrator パスワードは ADMIN です。小文字と大文字を切り替えるには仮想キーボードの Shift キーを使用します。

3. [Login  $( \mu \sigma / \gamma \nu )$ ] をタッチします。 これで Administrator アカウントに  $\mu \sigma / \gamma \nu \sigma$  ことができます。

- 注意 機器への無断アクセスを防ぐために、デフォルトのパスワードから新しいパスワードに変更して ください。パスワードの変更については、28 ページの「ユーザーアカウントを編集するには」を参 照してください。

#### 登録ユーザーとしてログインするには

- 1. ホーム画面で[Login (ログイン)]をタッチします。
- 2. Username ドロップダウンリストでユーザー名を選択し、パスワードを入力します。

#### ユーザーアカウントの管理

3. [Login (ログイン)]をタッチします。

Administrator は必要に応じてユーザーパスワードのリセットをすることができます。 28 ページの「ユーザーアカウントを編集するには」を参照してください。

#### ユーザーを変更するには

ログインするユーザーを変更するためにログアウトする必要はありません。

- 1. ホーム画面で[Login (ログイン)]をタッチします。
- 2. Username ドロップダウンリストメニューで変更したいユーザー名を選択します。

3.登録ユーザーとして、またはAdministratorとしてログインする場合にはパスワードを入力します。 ゲストとしてログインする場合には Password のテキストボックスを空白のままにします。

4. [Login (ログイン)]をタッチします。

確認のメッセージが表示されます。

### ユーザーアカウントの管理

管理者はホーム画面の Settings メニューを使用して、新規ユーザーの設定、ユーザーアカウント情報の編集、およびユーザーアカウントの削除を行うことができます。

新規ユーザーアカウントを追加するには

- 1. Administrator 権限でログインします。
- 2. ホーム画面で[Settings (設定)]をタッチします。

8	User Management	
Name	Access Level	Add
Administrator	Administrator	Add
Guest	Guest	
user123	Administrator	Edit
user456	User	
		Remove
		Cancel

[User Management (ユーザー管理)]をタッチします。
 User Management (ユーザー管理) 画面が開いて、使用できるユーザー名と対応するアクセスレベルが一覧で表示されます。

8	Add a New User
User Name Password Confirm Password	Access Administrator  Level

- 4. [Add (追加)]をタッチし、[Add a New User (新規ユーザーを追加)]画面を開きます。
- 5. [User Name (ユーザー名)]フィールドでユーザー名を入力します。
- 6. [Password (パスワード)]と[Confirm Password (パスワードの確認)]フィールドに新規アカウン ト用のパスワードを入力します。
- 7. [Account Level (アカウントレベル)]ドロップダウンリストでアカウントレベル(user または Administrator)を選択します。
   [User (ユーザー)]を選択すると登録ユーザーアカウントを作成します。
   [Administrator (管理者)]を選択すると、管理者権限を持つアカウントを作成します。

8、[OK]をタッチしてユーザーアカウントを保存します。

セットアップが成功すると確認のメッセージが表示され、登録したユーザー名が[User Management (ユーザー管理)]リストに追加されます。

#### ユーザーアカウントを編集するには

- 1. Administrator 権限でログインします。
- 2. ホーム画面で[Settings (設定)]をタッチします。
- [User Management (ユーザー管理)]をタッチします。
   User Management (ユーザー管理) 画面が開いて、使用できるユーザー名と対応するアクセス レベルが一覧で表示されます。
- 4. 編集したいユーザーアカウントを選択し、[Edit (編集)]をタッチすると[Edit User (ユーザーの編集)] 画面が表示されます。

88	Ed	it User	
User Name	user456	Access User	~
Password	****		
Password			

- 5. [Edit User (ユーザーの編集)]画面では、User レベルのアカウントのパスワードとアクセ スレベルを変更することができます。管理者とゲストのアカウントは編集することができません。
- [OK]をタッチして変更を保存します。
   ユーザーデータが更新され、確認メッセージが表示されます。

#### ネットワーク設定とリモートアクセス

#### ユーザーアカウントを削除するには

Administrator (管理者) および Guest (ゲスト) アカウントは削除することができません。

- 1. Administrator 権限でログインします。
- 2. ホーム画面で[Settings (設定)]をタッチします。
- [User Management (ユーザー管理)]をタッチします。
   User Management (ユーザー管理) 画面が開いて、使用できるユーザー名と対応するアクセス レベルが一覧で表示されます。
- 4. 削除するユーザー名を選択します。
- [Remove(削除)]をタッチすると選択したユーザーが削除されます。
   そのアカウントのユーザー権限が削除されます。また、そのユーザーに関連したすべてのプロ グラムと GLP レポートが削除されます。

# ネットワーク設定とリモートアクセス

#### 装置の IP アドレスの確認・編集

- 1・ホーム画面から[Settings (設定)]をタッチします。
- 2. 設定画面で[LAN Setting (LAN 設定)]をタッチします。装置がネットワークに接続されていれ ば LAN Settings 画面に IP アドレス、サブネットマスク、デフォルトゲートウェイ、DHCP サーバ ーが表示されます。

IP アドレスの取得には自動取得と手動設定の 2 通りの方法があります。

- [Obtain an IP Address automatically (IP アドレスの自動取得)]オプションを選択した場合、
   ネットワーク設定を入力するフィールドは使用しません。
- ・[Use Manual Configuration (手動設定を使用)]オプションを選択した場合、ネットワーク設定 のフィールドが入力可能になるので、必要な情報を入力します。ネットワーク設定を変更したら [OK]をタッチして保存します。

ログインすると、PCからウェブベースで SureCycler 8800 にアクセスして装置をコントロールする ことができます。PC 上の画面表示と、装置タッチスクリーン上の画面表示は若干異なります。

# 別の装置からのアクセス

- 1. ホーム画面で[Network Control (ネットワークコントロール)]をタッチします。
- 2、画面で6つの[Touch to Connect (タッチして接続)]の四角形のどれか1つをタッチします。

Touch to connect	Touch to connect	Touch to connect
Touch to connect	Touch to connect	Touch to connect

新しい画面が開きます。

- 3. 接続する装置の IP アドレスまたはサイクラー名を[IP Address/Cycler (IP アドレス/サイク ラー) フィールドに入力します。
- 4. [Connect (接続)]をタッチします。再び[Network Control (ネットワークコントロール)] 画面が表示され、接続する装置のネットワーク情報とステータスが四角形の中に表示されます。
- 5. 接続する装置の四角形をタッチします。その装置へのログインを求めるメッセージが表示されます。

🔨 Log In	The targeted machine is being used by Administrator . You can login to take over the access.
User Name Log In	

6. ユーザー名とパスワードを入力してリモートでログインします。この状態で、リモートで接続 した装置をコントロールすることが可能です。

# PC から装置へのアクセス

1. PC でインターネットブラウザ (Microsoft Internet Explorer が推奨)を開きます。

2. URL のフィールドに接続したい装置の IP アドレスを入力します。SureCycler 8800 ソフトウェ アが PC 上で表示されます。

🔨 Log In	The targeted machine is being used by Guest . You can login to take over the access.
User Name Log In	

3. ユーザー名とパスワードを入力してログインします。

装置本体の設定

# 装置本体の設定

### ホットトップの設定

サーマルサイクラーのホットトップのオン・オフの設定をするには

- 1. ホーム画面で[Settings (設定)]をタッチします。
- 2. [Hot Top Settings (ホットトップの設定)]をタッチします。
- 3. Hot Top Settings のスクリーンでホットトップ使用の有無を選択します。
  - ・[Yes]を選択するとサーマルサイクラーの運転中にホットトップを使用します。
  - ・[No]を選択すると、ホットトップはオフになります。
- 4. [OK]をタッチして選択を確認します。

ホットトップをオフにした場合、PCR 反応、インキュベーション反応のいずれの実行中にもホット トップは加熱されません。ダッシュボード上に、ホットトップ無効のアイコンが表示されます。 ホットトップの設定はユーザーレベルにかかわらず変更可能です。



# サーマルサイクラーの名前

サーマルサイクラーの名前は Administrator が変更することができます。

- 1. ホーム画面から[Setting (設定)]をタッチします。
- 2. [System Settings (システム設定)]をタッチします。
- 3. [Change (変更)]をタッチすると新しい画面が表示されて新しい名前を入力できるようにしま す。
- 4. [OK]をタッチして変更を保存します。

### 日付と時刻

日付と時刻の設定は Administrator が変更することができます。

- 1. ホーム画面から[Setting (設定)]をタッチします。
- 2. [System Settings (システム設定)]をタッチします。

3. 日付を修正したい場合は[System Date (システムの日付)]にあるカレンダーを使用して正し い日付を選択します。

System Date								
<< July 2010 >>								
Sun	Mon	Tue	Wed	Thu	Fri	Sat		
27	28	29	30	1	2	3		
4	5	6	7	8	9	10		
11	12	13	14	15	16	17		
18	19	20	21	22	23	24		
25	26	27	28	29	30	31		
1	2	3	4	5	6	7		
Today	y : 7/3	0/201	D	. U	2			

4 ・時刻を変更するには[System Time (システムの時刻)]で上向き/下向きの矢印をタッチして 分・秒を調整します。



5. [Apply (適用)]をタッチして変更を確認します。



4. PCR プログラム

新規 PCR プログラムの作成	36
手動でプログラムを作成するには	36
PCR ウィザードを使用して PCR プログラムを作成するには	39
384 ウェル用に PCR プログラムを作成する際のヒント	41
コマンドの説明	42
プリインストールされたプログラムの使用	51
PCR プログラムの編集	53
PCR プログラムの実行	55
インキュベーションの実行	56
レポートの表示	57
トラブルシューティング	58

この章では、PCR プログラムの作成、実行、インキュベーションの実行、レポートの表示などについて説明しています

# 新規 PCR プログラムの作成

新規 PCR を作成するには2つの方法があります。リストから必要なコマンドを選んでカスタマイズ したプログラムを手動で作成する方法と、ウィザードを使用してプログラムを半自動的に作成する 方法となります。以下のセクションではこれら新規プログラムの作成方法について説明します。

# 手動でプログラムを作成するには

新規 PCR プログラムは Programming screen で作成することができます。左側のコマンドリストか ら必要なステップを選び、プログラムを組み立てていきます。右側の Program Window には選択・ 追加されたコマンドが実行する順に並びます。図1 は、Hot Start (ホットスタート) コマンドを プログラムに追加した際の画面の例を示しています。

図1 SureCycler 8800 ソフトウェアのプログラミング画面



以下のステップに従い、新規 PCR プログラムを作成します。

1・ホーム画面で[New Program (新規プログラム)]をタッチすると、プログラム画面が開きます。

- 2. 画面左側の Command List (コマンドリスト)には、利用可能なコマンドが一覧表示されます。 新規 PCR プログラムに使用するコマンドと、その使用順序を決定します。コマンドの詳細につ いては、42 ページの「コマンドの説明」を参照してください。
- 3. 以下の2つの方法のいずれかを用いて Command List (コマンドリスト)から Program Window (プログラム画面) に一番目のコマンドを追加します。

コマンドをタッチしてから、矢印マーク(下図)をタッチして転送



・コマンドをダブルタッチ

いずれの方法でも新規画面が開き、温度や時間などのパラメータを入力することができます。

- 4. 表示に従い、追加したコマンドに必要な設定を入力します。
- 5. [OK]をタッチして設定を確認するとプログラミング画面に戻ります。

6. ステップ3からステップ5を繰り返し、新規 PCR プログラムに必要なコマンドを追加していき ます。Program Window のコマンド順序は、上向きと下向きの矢印ボタンを使用することで変更でき ます。



- 7. PCR プログラムについての情報を入力するには(オプション)
- a. [Program Information (プログラムの情報)]ボタン(下図)をタッチします。



b. キーボードのアイコンをタッチして、タッチスクリーンの仮想キーボードを表示します。



- c. プログラムの説明を入力し、終了したら [OK]をタッチしてプログラミング画面に戻ります。
- 8. PCR プログラムを実行する前にファイルを保存します。
- a. [Save (保存)]ボタンをタッチし、[Save Program (プログラムの保存)]画面を開きます。



 b. キーボードのアイコンをタッチしてタッチスクリーンの仮想キーボードを立ち上げ、[Program Name (プログラム名)]フィールドにプログラム名を入力します。名前の入力が終わったら、 アイコンをもう一度タッチして仮想キーボードを閉じます。



- c. 画面の下部にあるディレクトリを使用して、プログラムの保存先フォルダを選択します。新規 フォルダを作成したい場合は[Create Folder (フォルダ作成)]ボタンをタッチします。
- d. [Save to User Folder (ユーザーフォルダに保存)]フィールドに保存したいフォルダ名が正し く入っていることを確認し、[Save]をタッチしてプログラムを保存します。
- e. ファイルが正常に保存されると、ダイアログが表示されるので[OK]をタッチしてプログラミン グ画面に戻ります。

- 9. プログラムをスタートするには
- a. プログラミング画面で[Start]をタッチします。
- b. 表示にしたがって反応溶液量をマイクロリットル単位で入力し、[OK]をタッチします。

注 PCR プログラムの編集は、53 ページの「PCR プログラムの編集」を参照してください。

### PCR ウィザードを使用して PCR プログラムを作成するには

新規 PCR プログラムを作成する際に、SureCycler に搭載の PCR ウィザード機能を使用することもで きます。ウィザードを用いると、5 ステップで簡単に PCR プログラムを作成することができます。

#### PCR ウィザードの開始

・ホーム画面で[PCR Wizard (PCR ウィザード)]ボタンをタッチします。

PCR ウィザードでの5ステップ

#### 1. DNA ポリメラーゼの選択

・リストから使用する DNA ポリメラーゼを選択し、[Next]ボタンをタッチします。選択可能な DNA ポリメラーゼは以下のとおりです。

- ・Paq5000 (ホットスタートまたはスタンダードバージョン)
- PfuUltra II
- Herculse II
- ・*PfuTurbo*(ホットスタートまたはスタンダードバージョン)
- Taq

ウィザードのアルゴリズムにより、選択した酵素にあわせたガイドラインに基づいて各ステップで パラメータが決定されます。(それぞれの酵素の推奨サーマルサイクル条件については、酵素に付 属のマニュアルを参照してください。)各ステップでのパラメータは後から変更することも可能で す。詳細については 53 ページの「PCR プログラムの編集」を参照してください。

#### 4. PCR プログラム

PCR ウィザードを使用して PCR プログラムを作成するには

#### 2. プライマーの情報

・プライマーの情報を以下の2つの方法のどちらかで入力し、終了したら[Next]ボタンをタッチします。

・プライマー配列を直接入力するか、または USB メモリから配列を読み込みます。

・配列情報を入れない場合はプライマーの Tm 値を入力します。

ここで入力したプライマー配列情報は、PCR プログラムのアニーリング温度を設定するために使われます。アニーリング温度は2つのプライマーの Tm 値の平均より 5℃低く設定されます。プライマーの Tm 値ではなく配列を入力した場合には、それぞれのプライマーの Tm 値は以下の式にしたがって計算されます。

 $Tm = 69.3 + (0.41 \times GC\%) - (650/N)$ 

GC%はプライマー配列中のGC含量、Nはプライマー配列の塩基数を表します。

#### 3. PCR 産物の長さ

・予想されるPCR産物の長さを入力し、その単位としてbpまたはkbを選択します。終了したら[Next] ボタンをタッチします。

ここで入力した値は、PCR プログラム中で伸張反応の持続時間を設定するために使われます。

#### 4. DNA ソースの情報

・テンプレート DNA のソースをリストから選択し、[Next]をタッチします。ソースの選択肢は以下 のとおりです。

- cDNA
- ・ゲノム DNA
- ・ベクターDNA

DNA ソースの情報は、ホットスタート/初期変性の持続時間、伸張反応の持続時間を設定するために 使われます。

- 5. プログラムの調整
- 最後にプログラムの調整を行います。

・GC%は 70%以上ですか? テンプレート DNA が 70%以上の GC 塩基対を含む場合にこの調整を行います。テンプレート の Tm 値が高くなる分が補正されます。

・アニーリングのステップにグラジエントを設定しますか?

PCR プログラムのアニーリングのステップにグラジエントを設定したい場合に使用します。 ウィザードはプライマーの Tm 値に基づいて適切なグラジエント範囲を設定します。

ウィザードに従ってここまでの5ステップを終了すると、入力した情報に基づいた PCR プログラムが作成され、プログラミング画面に表示されます。

この時点で、手動でプログラムを作成した場合と同様に、各ステップのパラメータを変更したり、ステ ップを削除したりすることも可能です。詳しくは 53 ページの「プログラムのパラメータを編集するに」 を参照してください。

各ステップのパラメータを確認・編集するには、プログラムウィンドウでステップをダブルタッチしま す。新しい画面が開き、そのコマンドについての温度・時間などの設定が表示されます。各コマンドの 詳細については 42 ページの「コマンドの説明」を参照してください。

### 384 ウェル用に PCR プログラムを作成する際のヒント

384 ウェルプレートで PCR を行う場合、酵素によっては PCR サイクル中の熱変性ステップの持続時間を 短くすることにより、PCR 産物の収量が改善される場合もあります。プロトコルにより、3~5 秒の熱変 性時間が適している場合もあります。

PCR ウィザードを用いて PCR プログラムを作成する場合、プログラムが作成されてから熱変性の持続時間を短く変更する必要があります。詳しくは、53 ページの「パラメータの編集方法」を参照してください。

# コマンドの説明

以下のプログラムステップが PCR プログラムを作成する際に利用することができます。これらプロ グラムステップは、プログラミング画面左側のコマンドリストから選択できるようになっています。 それぞれのステップについて、設定が必要なパラメータが解説されています。

### Hot Start (ホットスタート)

ホットスタートプログラムステップは、ホットスタート仕様の DNA ポリメラーゼを使用する際の PCR プログラムの最初に設定される熱変性ステップです。この熱変性ステップで、DNA ポリメラー ゼに施された化学修飾をはずしたり、DNA ポリメラーゼに結合した抗体を変性させることにより、 DNA ポリメラーゼが活性化されます。ホットスタートは、ホットスタート仕様ではない DNA ポリメ ラーゼを使用する際にも、テンプレート DNA の初期変性のためとして使用される場合もあります。

ホットスタートのステップを追加する際には、温度と保持時間を設定する必要があります。プログ ラムをランさせた場合は、ホットスタートステップの直後に次のコマンドが実行されます。

🚱 Hot Start						
Hot Start Temperature	∞					
Hold For	00 : 00 min:sec					

Hot Start Temperature(ホットスタート温度):ホットスタートステップの温度を入力します。 Hold for (保持時間):ホットスタートステップの保持時間を入力します。適切に設定するため にDNA ポリメラーゼの製造元から提供される情報を参照してください。

# Cycle Wizard (サイクルウィザード)

サイクルウィザードコマンドは、複数回繰り返すサイクルを設定する際に使用します。サイクルウ ィザードを使うと、Start Cycle (サイクル開始)、End Cycle (サイクル終了) コマンドを手動 で追加する必要がなくなり、これらのコマンドがウィザードによって自動的に追加されます。

*	Cycle Wi	zard
Denaturation Step	95 °C	00:01:00 hour:min:sec
Annealing Step	62 ℃	00:00:30 hour:min:sec
✓ Use 3rd step Elongation Step	° 72℃	00:01:00 hour:min:sec
Cycles	40	

Denaturation Step (熱変性ステップ): 左側のフィールドに温度を℃単位で入力し、右側のフィールドに保持時間を時間:分: 秒の形式で入力します。

Annealing Step (アニーリングステップ): 左側のフィールドにアニーリング温度を℃単位で入力し、右側のフィールドに保持時間を時間:分:秒の形式で入力します。

Elongation Step (伸張ステップ): PCR プロトコルでアニーリングステップと伸張ステップを別々 にとる 3-step を採用する場合、Use 3-step 横のボックスにチェックを入れます。その上で、左側 のフィールドに温度を℃単位で入力し、右側のフィールドに保持時間を時間:分:秒の形式で入力 します。

Cycles (サイクル数): サイクル数を入力します。

# Start Cycle (サイクル開始)

このプログラムステップは PCR のサイクルを開始するために使用します。複数回繰り返したいステップの前にこのステップを置きます。

Start Cycle					
Number Of Cycles	40 cycles				
Cycle Name (Optional)	3-step with 62 degree annealing				

Number of Cycles (サイクル数): サイクルの繰り返し回数を入力します。

**Cycle Name**(サイクル名、オプション):必要に応じて、サイクルにも名前をつけることができます。

注

有効なプログラムを作成するためには、サイクルに含まれるステップは Start Cycle と

End Cycleのコマンドの間に入れる必要があります。

# Temperature Step (温度ステップ)

このプログラムステップでは、サーマルブロックを指定した温度で指定した時間保持します。さらに、温度や保持時間をサイクルごとに増加あるいは減少させることも可能です。また、サーマルブロックが設定温度に達する際の Ramp rate (ランプ・レート)を2段階の設定から選択することができます。

Temperature Step					
Processing Temperature	[55℃				
Hold For	00 : 0 : 30 hour:min:sec				
	Infinite				
Temperature Ramp	○ Standard				
Step Type	Annealing ~				
Increments per Cycle (Option	nal)				
Temperature Increment	C/cycle +/-				
Time Increment	5 seconds/cycle				

Processing Temperature (処理温度): ステップの温度を℃単位で入力します。

Hold for (保持時間): ステップの保持時間を時間:分:秒の形式で入力します。

**Temperature rate** (温度変化のレート): サーマルブロックの Ramp rate を選択します。

Fast ramp rate (高速ランプ・レート) は本機のデフォルトの Ramp rate で、大部分の PCR プロト コルに適したレートです。Standard ramp rate (スタンダード・ランプ・レート) はオプションと して用意されており、Standard ramp rate にすることで増幅が改善される場合もあります。Fast ramp rate を使用することで、アジレントの Herculase II DNA ポリメラーゼ、*PfuU/tra* II DNA Polymerase などの酵素のパフォーマンスが向上します。

**Step Type**(ステップのタイプ):ドロップダウンリストからステップのタイプを選択します。 Denaturation (熱変性)、Annealing (アニーリング)、Elongation (伸長)の3タイプから選択可 能です。 注 Increment per Cycle (サイクルごとの増分) フィールドは、サイクル中の温度ステップで 温度、時間をサイクルごとに増減させる設定のために使用します。この増減設定はオプション ですので不要な場合は入力せずにフィールドを空白のままにしてください。

Temperature Increment (温度増分): サイクルごとの温度の増分を入力します。+/-ボタンをタッチして値の正負を切り替えます。負の値(たとえば-2℃)はサイクルごとに温度を下げることを意味します。

**Time Increment** (時間増分): サイクルごとの Hold for (保持時間)の増分を入力します。この フィールドの単位は秒/サイクルです。

# End Cycle (サイクル終了)

PCR サイクルを終了させるためにこのプログラムステップを挿入します。複数回繰り返したい一連のステップのあとにこのステップを置きます。

**注** 有効なプログラムを作成するためには、サイクルに含まれるステップは Start Cycle (サイク

ル開始)とEnd Cycle(サイクル終了)のコマンドの間に入れる必要があります。

# Gradient Step (グラジエントステップ)

Gradient Step (グラジエントステップ)は Temperature Step (温度ステップ) に類似しています が、Temperature step ではブロック全体を同一の温度で保持するのに対して Gradient Step ではサ ーマルブロックに温度グラジエントをかけて指定の時間、保持します。

Gradient Step は特定のプライマー・テンプレートに対してアニーリング温度を最適化する場合に 便利です。

#### 4. PCR プログラム

Gradient Step (グラジエントステップ)

ランの実行中、Left Side Temperature Gradient(グラジエントの左端温度)と Right Side Temperature Gradient(グラジエントの右端温度)に入力された値に基づいてサーマルブロックに は温度グラジエントが左右方向にかかります。グラジエントの最低温度と最高温度の差が 30℃を超 える設定はできません。

Gradient Step													
Left Side Temperature Gradient	58	1	2	m	4	Li	و	7		e	9	:	B
Right Side Temperature Gradient	66	28.0 °C	58.7 °C	Do 2.62	60.2 °C	Do 6.03	61.6 °C	62.4 °C	63.1 °C	63.8 °C	64.5 °C	65.3 °C	66.0 °C
Hold For	00:30	m	in:	sec	2								
Convert to Temperature Step (optional)  After optimization, convert this gradient step of the program to a standard temperature step using the temperature from column: 1													

Left Side Temperature Gradient (グラジエントの左端温度): サーマルブロックの左端 (列番号 1) に設定したい温度を入力します。

**Right Side Temperature Gradient**(グラジエントの右端温度): サーマルブロックの右端(列番号 12) に設定したい温度を入力します。

Hold for (保持時間): グラジエントステップの保持時間を分: 秒の形式で入力します。

Convert to Temperature Step (optional) (温度ステップへの変換 オプション): 最適なアニ ーリング温度を決定したら、グラジエントステップから通常の温度ステップへと簡単に変換するこ とができます。もっとも増幅結果のよかった列が何列目だったかをドロップダウンリストから選択 すると、サーマルブロック全体がその列の温度設定となる温度ステップへと自動変換されます。

# Touchdown (タッチダウン)

タッチダウン PCR はプライマーとテンプレートの非特異的な結合を減らすための応用法です。タッ チダウン PCR では初期のサイクルでアニーリング温度を比較的高く設定し、特異的な増幅を促進さ せます。それ以降のサイクルではアニーリング温度を下げることでより増幅しやすい条件にします。

タッチダウンプログラムステップは、タッチダウン PCR を簡単にプログラミングできる1ステップ のコマンドです。アニーリング温度の最高値(サイクルスタート時のアニーリング温度)と、最低 値(最終サイクルのアニーリング温度)を入力します。ソフトウェアは各サイクルで均等にアニー リング温度が下がっていくように計算し、ランを実行します。

Touchdown						
Denaturation Temperature	95 °C					
Hold For	00 : 30 min:sec					
Max Annealing Temperature	65 ℃					
Min Annealing Temperature	60 °C					
Hold For	00 : 30 min:sec					
Elongation Temperature	[72℃					
Hold For	01 : 00 min:sec					
Cycles	40					

**Denaturation Temperature and Hold Time** (熱変性温度・保持時間): 熱変性温度を Denaturation Temparature フィールドに入力し、保持時間をその下の Hold For フィールドに入力します。

Max Annealing Temperature (最高アニーリング温度):アニーリング温度の最高値を入力します。 PCR の最初のサイクルのアニーリング温度にこの値が使用されます。

Min Annealing Temperature (最低アニーリング温度):アニーリング温度の最低値を入力します。 PCR の最後のサイクルのアニーリング温度にこの値が使用されます。

#### 注

アニーリング温度の最高値と最低値の差は通常 5~10°C程度で、Tm 値が高いほうのプライマー

の Tm 値よりも 2℃高い温度を最高温度として反応を開始します。たとえば、プライマー1の Tm 値が 60℃、プライマー2 の Tm 値が 54℃の場合はアニーリング温度の最高値は 62℃とし、 最低値は 52℃あたりに設定します。

Hold For (保持時間): アニーリングステップの保持時間を入力します。 Elongation Temperature and Hold Time (伸長反応の温度・保持時間): タッチダウンサイクル中 に使用する伸長反応の時間および保持時間を入力します。 Cycles (サイクル数): タッチダウンステップのサイクル数を入力します。 Strage / Pause Step (保存/一時停止ステップ)

# Storage/Pause Step (保存/一時停止 ステップ)

Storage/Pause(保存/一時停止)ステップでは、指定した時間または無期限にサーマルブロックを 4~12℃の範囲で冷却することができます。

Storage / Pause Step					
Hold Temperature 4 °C					
Hold For 02 : 00 : 00 hour:min:sec					
O Infinite					

Hold Temperature (保持温度):保持温度を4~12℃の範囲で入力します。

**Hold for** (保持時間):時間を指定する場合は Hold for のラジオボタンを選択し、保持時間を時間:分:秒の単位で入力します。

**Infinite** (無期限):時間を指定せずに無期限で保持したい場合には、Infiniteのラジオボタン を選択します。 4. PCR プログラム プリインストールされたプログラムの使用

# プリインストールされたプログラムの使用

SureCycler 8800 には 96 ウェルブロックで使用可能な PCR プログラムがあらかじめインストールされています。これらのプログラムはアジレントの PCR 酵素や変異導入キットにあわせてデザインされています。

それぞれのプログラムは、特定の長さのターゲットを一般的なアニーリング条件で増幅する場合に 合わせて設定されています。ターゲットの長さやプライマーのTm値によってはアニーリング温度、 伸長時間を調節する必要があります。プリインストールされたプログラムはどれも別のファイル名 で保存し、編集することができます。

プリインストールされているプログラムを表5に示します。

表5 プリインストールされた 96 ウェル SureCycler 8800 用 PCR プログラム

PCR Program File Name	PCR Enzyme or Kit	Denaturation	Annealing	Elongation	
Herculase II Fusion 1kb target.pcr	Herculase II Fusion DNA Polymerase	culase II Fusion DNA Polymerase 95°C for 20 seconds			
Paq5000 HS MM 1kb target.pcr	Paq5000 Hotstart Master Mix	95°C for 20 seconds	55°C for 20 seconds	72°C for 30 seconds	
PfuTurbo HS MM 1kb target.pcr	PfuTurbo Hotstart Master Mix	95°C for 30 seconds	55°C for 30 seconds	72°C for 1 minute	
PfuUltra II HS MM 1kb target.pcr PfuUltra II Hotstart Master I		95°C for	55°C for	72°C for	
		20 seconds	20 seconds	15 seconds	
QuikChange II 4 kb target.pcr	QuikChange II Site-Directed	95°C for	55°C for	68°C for	
	Mutagenesis Kit	30 seconds	1 minute	4 minutes	
QuikChange Lightning 4kb target.pcr	QuikChange Lightning Site-Directed	95°C for	60°C for	68°C for	
	Mutagenesis Kit	20 seconds	10 seconds	2 minutes	
QuikChange Lighting Multi 4 kb	QuikChange Lightning Multi	95°C for	55°C for	65°C for	
target.pcr	Site-Directed Mutagenesis Kit	20 seconds	30 seconds	2 minutes	
QuikChange Multi 4kb target.pcr	QuikChange Multi Site-Directed	95°C for	55°C for	65°C for	
	Mutagenesis Kit	1 minute	1 minute	8 minutes	

# プリインストールされたプログラムの利用

- 1. ホーム画面で[Program (プログラム)]をタッチします。
- 2. 画面左側のブラウザで Guest/Agilent フォルダに移動します。 プリインストールされたプログ ラムはこのフォルダに保存されています。
- (使用するプログラムをタッチして選択します。選択したプログラムを実行するには[Run(実行)]
   をタッチします。選択したプログラムのパラメータを確認するには[Open(開く)]をタッチします。

### プリインストールされたプログラムの編集

- 1. ホーム画面で[Program (プログラム)]をタッチします。
- 2. 画面左側のブラウザで Guest/Agilent フォルダに移動します。
- 3. 編集するプログラムを選択し[Open (開く)]をタッチします。プログラムがプログラミング画 面で開きます。
- 4. プログラミング画面で[Save (保存)]をタッチし、ファイルを別名で保存します。
- 5. プログラムを編集します。編集方法は 53 ページの「PCR プログラムの編集」を参照してください。

# PCR プログラムの編集

手動で作成したプログラム、ウィザードを使用して作成したプログラムはいずれもプログラミング 画面の機能を使って編集することができます。プリインストールされたプログラムもいったん別名 で保存すれば同様に編集することができます。

PCR プログラムを編集した後にプログラムは同じファイル名のまま、あるいは別のファイル名で保 存することができます。USB メモリに PCR プログラムをバックアップしたり、USB メモリを用いて 別の SureCycler 8800 にプログラムを転送することもできます。

### プログラムのパラメータを編集するには

- ホーム画面で[Programs (プログラム)]をタッチします。
   ユーザーレベル(ゲストまたは登録ユーザー)に応じて、特定のユーザーディレクトリまたはゲ ストディレクトリが表示され、既存の PCR プログラムがすべて表示されます。
- 編集するプログラムを選択して[Open(開く)]をタッチします。
   プログラミング画面の Program Window(プログラムウィンドウ)パネルにプログラムが表示されます。
- 3. 編集したいステップをダブルタッチすると、プログラムステップの編集画面が開きます。
- 4. 必要に応じてパラメータを編集し、[OK]をタッチするとプログラミング画面に戻ります。
- 5. ステップ3とステップ4を繰り返して必要な設定変更を行います。
- 6. プログラミング画面で[Save(保存)]をタッチします。同じプログラムのまま変更を保存する か、新しいプログラム名をつけて保存することができます。
- 7. ここで[Start (開始)]をタッチすれば PCR プログラムが開始されます。

### PCR ステップをコピーするには

- 1. プログラミング画面に移動します。
- 2. Program Window (プログラムウィンドウ) でコピーしたいステップを選択します。
- 3. コピーボタン(下記)をタッチします。



### PCR ステップを追加または削除するには

1. プログラミング画面に移動します。

新規ステップを追加するには、画面左側の[Command List(コマンドリスト)] でコマンドをタッチ して、右向きの矢印ボタンをタッチします。表示に従い、追加したコマンドに必要な設定を入力し ます。各コマンドの詳細は 42 ページの「コマンドの説明」を参照してください。

プログラムからステップを削除するには、Program Window(プログラムウィンドウ)内で削除した いステップを選択し、[Delete(削除)](下記)をタッチします。



プログラムの変更を保存するには、[Save(保存)]をタッチします。同じプログラムのまま変更を 保存するか、新しいプログラム名をつけて保存することができます。

### PCR プログラムを削除するには

- ホーム画面で[Programs (プログラム)]をタッチします。ログインしているユーザーレベル に応じて (ゲストまたは登録ユーザー)、特定のユーザーフォルダあるいはゲストフォルダに既 存の PCR プログラムが表示されます。
- 2. 削除したいプログラムを選択します。
- 3. [Delete(削除)]をタッチしてプログラムをフォルダから削除します。

# PCR プログラムの実行

# PCR プログラムを実行するには

ホーム画面からプログラムを実行するには

- 1. [Programs (プログラム)]をタッチします。
- 2. 実行したいプログラムをタッチします。
- 3. [Run Selected (実行)]をタッチします。

4. ダイアログボックスに反応溶液のボリュームをマイクロリットル単位で入力し、[OK]をタッチ します。

#### プログラミング画面からプログラムを実行するには

1. [Start]をタッチします。

2. ダイアログボックスに反応溶液のボリュームをマイクロリットル単位で入力し、[OK]をタッチ します。

### 実行中のプログラムを中止または一時停止するには

ステータス画面から:

[Stop(停止)](以下に表示)をタッチし、プログラムの実行を中止します。



[Pause(一時停止)](以下に表示)をタッチし、プログラムを一時停止します。[Pause(一時停止)]はプログラムの実行中にのみ有効になります。一時停止すると、Pause ボタンは Play ボタン にかわります。プログラムを再開するには Play ボタンをタッチします。



#### **4. PCR プログラム**

インキュベーションの実行

# インキュベーションの実行

インキュベーション機能は指定した時間または無期限に、一定の温度を保ってインキュベーション を行う際に便利です。

# インキュベーションの設定と実行

- 1. ホーム画面で[Incubation]をタッチします。
- 2. [Incubation Temperature]フィールドでインキュベーション温度を入力します。
- 3. インキュベーションの保持時間を指定します。

・指定した時間、一定温度で保持する場合は[Hold for (保持時間)]を選択し、時間:分:秒の 形式で保持時間を入力します。

- ・無期限に保持する場合は、[Infinite(無期限)]を選択します。
- 4. [OK]をタッチしてインキュベーションを開始します。

# レポートの表示

PCR プログラムが実行されると、GLP レポートが作成されユーザーフォルダに保存されます。たと えば、ゲストアカウントで PCR プログラムを実行すると、GLP レポートはゲストフォルダに保存さ れます。GLP レポートのファイルは.glp という拡張子を持ちます。

### レポートを表示するには

- ホーム画面で[Reports (レポート)]をタッチします。
   レポートブラウザが起動し、レポートファイルのリストが表示されます。
- スクリーン下の矢印ボタンを使用してレポートのリストをスクロールすることができます。また、レポートを名前や日付でソートすることもできます。[Report Name]カラムのヘッダーをタッチするとレポートの名前でソートされ、[Date]カラムのヘッダーをタッチするとレポートの日付でソートされます。タッチスクリーン上でドラッグすることにより、カラムの幅を変えることができます。
- 3. 表示させたいレポートを選び、[Open]をタッチします。レポートは新しいスクリーンで表示さ れます。[Back]ボタンをタッチすると、レポートブラウザに戻ります。

### レポートを削除するには

レポートブラウザで削除するレポートを選択し、[Delete(削除)]をタッチします。

トラブルシューティング

# スタンダード・ランプレートを使って、少量で反応させた際に PCR 産物の収

#### 量が低い場合

- ✓ プレートをシーリングフィルムで閉じ、コンプレッション・マットを使用している場合、サイクルの中の熱変性の時間を短縮します。熱変性時間は 3~5 秒程度で適切な場合があります。 または、シーリングフィルムではなくキャップを使用します。
- ✓ 反応ボリュームを 50 uL に増やすと改善される場合があります。
- ✓ ファスト・ランプレートへ切り替えることで改善される場合があります。

### プログラムの実行中にブロック温度が表示されない場合

サーマルブロックユニットが正しくセットされていない可能性があります。

✓ 電源を抜いた上でサーマルブロックユニットが正しく取り付けられていることを確認し、装置 を再起動します。

### プログラムの実行が始まらない場合

サーマルブロックユニットが正しく取り付けられていない可能性があります。

✓ 電源を抜いた上でサーマルブロックユニットが正しく取り付けられていることを確認し、装置 を再起動します。

# チューブの内部に結露があった場合

PCR の終了時に結露するのは自然な現象であり、有害な効果はありません。

✓ チューブをスピンダウンして結露した水分を反応液中に戻した上で、その後の実験に用いてく ださい

ホットトップが無効になっていると水分が蒸発しやすくなります。

✓ ホットトップを有効にしてください。詳細は 32 ページの「ホットトップの設定」を参照して ください。

### シールしたマイクロプレートでサンプル液量が減少した場合

- ✓ マイクロプレート、シールの品質が十分に良いものであることを確認してください。
- ✓ PCR 中にシーリングフィルムがプレートに密着した状態が保たれるように、コンプレッション・マットを使用してください。

### メモリスティック、キーボードなどの USB デバイスが認識されない場合

✓ 装置を再起動してください。

#### 装置の電源が入らない場合

- ✓ 電源コードが正しく接続され、電力が供給されていることを確認してください。
- ✓ 装置後部の電源コード差込口の上に位置するヒューズが切れていないか確認してください。

# 画面が点灯せず、何も表示されない場合

- ✓ 電源コードが正しく接続され、電力が供給されていることを確認してください。
- ✓ 装置後部の電源コード差込口の上に位置するヒューズが切れていないか確認してください。

# 画面が点灯するものの、何も表示されない場合

- ✓ 装置を再起動してください
- ✓ テクニカルサポートに連絡してください。国別の連絡先は14ページに記載されています。



5. メンテナンスと ソフトウェアの更新

メンテナンス	62
クリーニング	62
ヒューズの交換	62
ソフトウェアの更新	63

この章では、機器のメンテナンスおよびソフトウェア更新のアップロードに関する方法について説明しています。

5. メンテナンスとソフトウェアの更新 メンテナンス

# メンテナンス

### クリーニング

SureCycler 8800 のメンテナンス方法を示します。

・クリーニング前に必ず本体の電源をオフにするか、電源を外してください。

・本体およびサーマルブロックのウェルの清掃には、水またはイソプロピルアルコールで布や綿棒などを用いて行ってください。

・その他の有機溶媒はSureCycler 8800の清掃には用いないでください。

・液体が本体内に入らないようにご注意ください。

# ヒューズの交換

**警告** 感電を防ぐため、ヒューズ交換を行う前に必ず電源コードの接続を外してください。

ヒューズコンパートメントは本体背面の主電源接続部の上にあります。ヒューズの規格が T10A、 250V であることをチェックしてください。

### ソフトウェアの更新

SureCycler 8800 ソフトウェアの新バージョンがリリースされると、アジレントのウェブサイトか らダウンロードし、ご利用の SureCycler 8800 にアップロードすることができます。

- 新しいバージョンのソフトウェアがご利用可能になった際、弊社よりお知らせを通知いたします。このお知らせには、ソフトウェアのファイルにアクセスするためのウェブアドレスが記載されています。ダウンロードしたファイルは USB メモリスティックに保存し SureCycler 8800本体にインストールすることが可能です。
- SureCycler 8800本体の Administrator アクセス権限を持つユーザーアカウントにログインします(操作方法については25ページの「管理者としてログインするには」をご参照ください)。
- 3. ホーム画面で[Settings (設定)]、[Software Update (ソフトウェアの更新)]をタッチしま す。
- 本体前面にある USB ポートのどちらかにソフトウェアを保存した USB メモリスティックを挿入 します。
- [Browse (ブラウズ)]ボタンをタッチしてブラウザを開きます。ブラウザでソフトウェアが保存されているフォルダまで移動します。適切なフォルダを選択して[OK]ボタンをタッチすると、 [Software Update (ソフトウェアの更新)]画面に戻ります。
- 6. [Update(更新)]をタッチします。選択したフォルダでソフトウェアファイルが認識されてい ない場合には[Update(更新)]ボタンを利用することができません。
- 本体へのファイルのダウンロードが開始されます。すべてのファイルのダウンロードが完了すると、自動的に再起動します。
- 8. 再起動が開始したら、USBメモリスティックを本体から取り外します。
- 9. 再起動が完了したら、[Software Update (ソフトウェアの更新)]画面に戻って新しいバージョンのソフトウェアが動作していることを確認してください。

# 本書では

本文書は Agilent SureCycler 8800 の設置方法、使用方法について説明 しています。

© Agilent Technologies, Inc. 2011

バージョンA, 2011 年 3 月



G8800-95000

