

# Applied Biosystems Veriti™ サーマルサイクラー

日本語ユーザーマニュアル



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。

This Veriti<sup>™</sup> Thermal Cycler is covered by patents owned by Applera Corporation. No right is conveyed expressly, by implication, or by estoppel to apparatus, reagents, kits, or methods such as the patented 5' nuclease methods. For further information contact the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

Applied Biosystems, BigDye and MicroAmp are registered trademarks and AB (Design), Applera, MicroAmp, Veriti and VeriFlex are trademarks of Applera Corporation or its subsidiaries in the US and/or certain other countries.

本誌に記載の社名および製品名は、弊社または各社の商標または登録商標です。

記述内容は改良のため予告なく変更される場合がありますので、ご了承下さい。

製品の仕様・外観は改良のため予告なしに変更される場合があります。

ご注意

この簡易ユーザーマニュアルは、機器付属のユーザーマニュアル(P/N 04375799)を簡略化 したものです。

詳細な説明や知的所有権等に関しては付属のユーザーガイドを必ずご確認ください。

© 2007 Applied Biosystems Japan Ltd,. All rights reserved.

目 次

1.	システムの概要	4
	1.1 Veriti <sup>™</sup> サーマルサイクラーについて	5
	装置について	5
	1.2 タッチスクリーンの利用	6
	文字の入力	7
	数字の入力	8
	1.3 メニューの概要	9
	Main Menu 画面	9
	Settings Menu 画面	10
	Admin Menu 画面	10
	Tools Menu 画面	11
	1.4 ランメソッドについて	11
	ランメソッド とは	11
	デフォルトで 保存されているランメソッド	11
	ランメソッドの 閲覧	12
	ランメソッドの 管理	13
	ランメソッドの コピー	14
	ランメソッドの 消去	15
	ランメソッドを デフォルト として設定	15
	ランメソッド 及び各設定の バックアップ	16
	ランメソッドや その他の設定の 回復・転送	17
	1.5 ランフォルダについて	18
	カスタムフォルダの消去	18
	カスタルフォルダの名称亦再	10
	77774777777777777777777777777777777777	19
2.	ランの実行	20
2.	ランの実行 2.1 消耗品の選択	19 20
2.	<b>ランの実行</b> 2.1 消耗品の選択 チューブの選択	19 20 21
2.	<ul> <li>ランの実行</li> <li>2.1 消耗品の選択</li></ul>	19 20 21 21 21
2.	<ul> <li>ランの実行</li> <li>2.1 消耗品の選択</li></ul>	19 20 21 21 21 22 26
2.	<ul> <li>ランの実行</li> <li>2.1 消耗品の選択</li></ul>	19 20 21 21 21 22 26 26
2.	<ul> <li>ランの実行</li> <li>2.1 消耗品の選択</li></ul>	19 20 21 21 22 26 26 26
2.	<ul> <li>ランの実行</li> <li>2.1 消耗品の選択</li></ul>	19 20 21 21 22 26 26 26 26
2.	<b>ランの実行</b> 2.1 消耗品の選択 チューブの選択 サンプルトレーとプレートの 組み合わせ 2.2 サンプルを装置にセットする Veriti サーマルサイクラー: 0.2mL ブロックの場合 キャップ付き チューブの セット 96-well プレートのセット キャップなし チューブの セット	19 20 21 21 22 26 26 26 26 26 26 27
2.	<b>ランの実行</b> 2.1 消耗品の選択 チューブの選択 サンプルトレーとプレートの 組み合わせ 2.2 サンプルを装置にセットする Veriti サーマルサイクラー: 0.2mL ブロックの場合 キャップ付き チューブの セット 96-well プレートのセット キャップなし チューブの セット Veriti Fast サーマルサイクラー: 0.1mL ブロックの場合	19 20 21 21 22 26 26 26 26 26 26 27 27
2.	<b>ランの実行</b> 2.1 消耗品の選択 チューブの選択 サンプルトレーとプレートの 組み合わせ 2.2 サンプルを装置にセットする Veriti サーマルサイクラー: 0.2mL ブロックの場合 キャップ付き チューブの セット 96-well プレートのセット キャップなし チューブの セット Veriti Fast サーマルサイクラー: 0.1mL ブロックの場合 キャップ付き チューブの セット	19 20 21 21 22 26 26 26 26 26 26 27 27 27
2.	<b>ランの実行</b> 2.1 消耗品の選択 チューブの選択 サンプルトレーとプレートの 組み合わせ 2.2 サンプルを装置にセットする Veriti サーマルサイクラー: 0.2mL ブロックの場合 キャップ付き チューブの セット 96-well プレートのセット キャップなし チューブの セット Veriti Fast サーマルサイクラー: 0.1mL ブロックの場合 キャップ付き チューブの セット 96-well デューブの セット	19 20 21 21 22 26 26 26 26 26 26 27 27 27
2.	ランの実行 2.1 消耗品の選択	19 20 21 21 22 26 26 26 26 26 26 27 27 27 27 27
2.	ウンの実行 2.1 消耗品の選択 チューブの選択 サンプルトレーとプレートの 組み合わせ 2.2 サンプルを装置にセットする Veriti サーマルサイクラー: 0.2mL ブロックの場合 キャップ付き チューブの セット 96-well プレートのセット キャップなし チューブの セット Veriti Fast サーマルサイクラー: 0.1mL ブロックの場合 キャップ付き チューブの セット Set サーマルサイクラー: 0.1mL ブロックの場合 キャップ付き チューブの セット Neriti Fast サーマルサイクラー: 0.1mL ブロックの場合 キャップ付き チューブの セット 96-well Fast プレートの セット 96-well Fast プレートの セット 96-well Fast プレートの セット 96-well Fast プレートの セット	19 20 21 21 22 26 26 26 26 26 26 27 27 27 27 27 27 28 28
2.	<b>ランの実行</b> 2.1 消耗品の選択 チューブの選択 サンプルトレーとプレートの 組み合わせ 2.2 サンプルを装置にセットする Veriti サーマルサイクラー: 0.2mL ブロックの場合 キャップ付き チューブの セット 96-well プレートのセット キャップなし チューブの セット Veriti Fast サーマルサイクラー: 0.1mL ブロックの場合 キャップ付き チューブの セット Second ディーブの セット シートの セット シートの セット 96-well Fast プレートの セット 96-well Fast プレートの セット 96-well Fast プレートの セット 2.3 新しいランメソッドの作成	19 20 21 21 22 26 26 26 26 26 26 26 27 27 27 27 27 27 27 27 27 27 28 29
2.	<b>ランの実行</b> 2.1 消耗品の選択         チューブの選択         サンプルトレーとプレートの 組み合わせ         2.2 サンプルを装置にセットする         Veriti サーマルサイクラー: 0.2mL ブロックの場合         キャップ付き チューブの セット         96-well プレートのセット         キャップなし チューブの セット         Veriti Fast サーマルサイクラー: 0.1mL ブロックの場合         キャップなし チューブの セット         Se-well Fast ブレートの セット         96-well Fast ブレートの セット         Veriti Fast サーマルサイクラー: 0.1mL ブロックの場合         キャップはき チューブの セット         Se-well Fast ブレートの セット         96-well Fast ブレートの セット         93 新しいランメソッドの作成         ランメソッドの新規作成	19 20 21 21 22 26 26 26 26 26 26 27 28 28 29 29 27 
2.	<b>ランの実行</b> 2.1 消耗品の選択 チューブの選択 サンプルトレーとプレートの 組み合わせ 2.2 サンプルを装置にセットする Veriti サーマルサイクラー: 0.2mL ブロックの場合 キャップ付き チューブの セット 96-well プレートのセット キャップなし チューブの セット Veriti Fast サーマルサイクラー: 0.1mL ブロックの場合 キャップはき チューブの セット 96-well デレートの セット 96-well Fast プレートの セット 96-well Fast プレートの セット 93 新しいランメソッドの作成 ランメソッドの新規作成 2.4 ランメソッドの編集	19 20 21 21 22 26 26 26 26 26 26 27 27 27 27 27 27 27 28 28 29 29 31
2.	ランの実行         2.1 消耗品の選択         チューブの選択         サンプルトレーとプレートの組み合わせ         2.2 サンプルを装置にセットする         Veriti サーマルサイクラー: 0.2mL ブロックの場合         キャップ付き チューブの セット         96-well プレートのセット         キャップなし チューブの セット         Veriti Fast サーマルサイクラー: 0.1mL ブロックの場合         キャップなし チューブの セット         96-well プレートの セット         Po-well Fast プレートの セット         Po-well Fast プレートの セット         Stube         シンプルトレーまたはプレートをサンプルブロック にセット         2.3 新しいランメソッドの作成         ランメソッドの編集         ステージの追加	19 20 21 21 22 26 26 26 26 26 26 27 27 27 27 27 27 27 28 28 28 29 31 31
2.	<b>ランの実行</b> 2.1 消耗品の選択 チューブの選択 サンプルトレーとプレートの 組み合わせ 2.2 サンプルを装置にセットする Veriti サーマルサイクラー: 0.2mL ブロックの場合 キャップ付き チューブの セット 96-well プレートのセット キャップなし チューブの セット Veriti Fast サーマルサイクラー: 0.1mL ブロックの場合 キャップ付き チューブの セット Veriti Fast サーマルサイクラー: 0.1mL ブロックの場合 キャップ付き チューブの セット S6-well Fast プレートの セット 96-well Fast プレートの セット 2.3 新しいランメソッドの作成 ランメソッドの新規作成. 2.4 ランメソッドの編集 ステージの追加 ステップの追加	19 20 21 21 22 26 26 26 26 26 26 27 27 27 27 27 27 27 27 28 29 31 31 32

	温度の移行速度の変更	33
	VeriFlex ブロックで 異なる温度プログラムを設定	
	AutoDelta ステップの設定	
	ポーズの追加	
	ステージまたは ステップの消去	
	2.5 ランの実行	
	ランを開始する前に	
	2.6 ランのモニタリング	
	Run Monitor 画面	
	ラン中の Run Monitor 画面における操作	41
	サイクル数の変更	
	ランのポーズ	
	ランの停止	
	Run Monitor ボタンを使う	
	2.7 ランレポートの閲覧と印刷	
	2.8 装置からサンプルを取り出す	
	2.9 サイクルパフォーマンステストの実行	
	必要なもの	
	操作手順	
3.	Veriti <sup>™</sup> サーマルサイクラーの初期設定	46
	日付と時間の設定	46
	セキュリティおよびユーザーアカウントの設定	
	セキュリティ設定の有効化	
	アクセスレベル	
	新規ユーザーの 設定	
	ユーザーの追加	
	ユーザーの 変更・編集	
	ユーザーの削除	
	その他のオプション設定	
	ヒートカバー及びサンプルブロックのアイドリング 温度の設定	
	ポーズまたは スタンバイの 持続時間	
	Main Menu 画面の Shortcut の設定	53

## 1. システムの概要

このセクションでは下記について解説しています。

- 1.1 Veriti<sup>™</sup>サーマルサイクラーについて
- 1.2 タッチスクリーンの利用
- 1.3 メニューの概要
- 1.4 ランメソッドについて
- 1.5 ランフォルダについて

## 1.1 Veriti<sup>™</sup>サーマルサイクラーについて

装置について Veriti<sup>™</sup>サーマルサイクラーには2つのモデルがあります。

- Veriti<sup>™</sup> 96-Well Fast サーマルサイクラー、0.1mL サンプルブロック (商品番号 Veriti Fast 100)
- Veriti<sup>™</sup> 96-Well サーマルサイクラー、0.2mL サンプルブロック (商品番号 Veriti 200)

Veriti<sup>™</sup>サーマルサイクラーには"VeriFlex<sup>™</sup> サンプルブロック"という機能 が搭載されています。この機能はサンプルブロック上の独立した6つのエリア 毎に任意の温度設定が可能で、複雑なプライマーセットのデザインや PCR 条件 の最適化検討に役立ちます。また、大型カラータッチ画面による操作と分かり やすいインターフェイスにより、快適な環境で使用できます。 Veriti<sup>™</sup>サーマルサイクラーは、ネットワークを通じて最大12台まで接続する ことができ、これにより、ある装置のランメソッドを他の装置へ転送したり、 一度に複数台の装置を同時にランスタートすることができます。

このガイドには、Veriti<sup>™</sup>サーマルサイクラーの概要および基本的な操作について解説しています。



アプライドバイオシステムズ Veriti<sup>™</sup> サーマルサイクラーの外観

## 1.2 タッチスクリーンの利用

タッチ画面を使用することで、Veriti™サーマルサイクラーを快適に操作する ことが可能です。

下表にタッチ画面上でよく使うボタンについて説明しています。

ボタン 機能 Main Menu 画面に戻ります。 1つ前の画面に戻ります。 Done Х 装置が運転中に現れ、運転中に生じたエラー を表示するステータスレポート画面を表示 します。 装置が運転中に現れ、運転状況を表示する Run Monitor 画面を表示します。 Stage 1 Time Remaining: Sample: 61.0 °C ランメソッドなどのリストの表示項目を上 下にスクロールします。 ¥ ランメソッドフォルダを表示します。 D 装置の USB ポートに USB ドライブが接続 されていることを示しています。 ÷ ヘルプメニューを表示します。

Veriti<sup>™</sup> サーマルサイクラーのタッチ画面上のボタン

**文字の入力** ユーザー名やパスワードなどの文字入力が必要な場合、図 1-2 に示す文字入力 画面が現れます。



Veriti™サーマルサイクラーの文字入力画面

操作	使用するキー		
小文字を入力する	直接文字ボタンで入力する。		
大文字を入力する	Caps ボタンを押した後、文字ボ           タンを入力する。もう一度 Caps           ボタンを押すと、小文字入力画           面に戻ります。		
記号や句読点を入力する	Symbols か         Symbols ボタンを押した           後、記号ボタンで入力す         る。再度 Symbols ボタン           る。再度 Symbols ボタン         を押すと、小文字入力画           面に戻ります。         1		
入力した文字を全て消去する	Clear		
最後に入力した文字を消去する 			
入力画面を終了する	Done		

### 数字の入力 数字の入力が必要な場合、下図に示す数字入力画面が現れます。



Veriti™サーマルサイクラーの数字入力画面

操作	使用するキー
数字を入力する	直接数字ボタンで入力します。 この装置は温度の入力など小数点を含 むフィールドでは、自動的に小数点を表 示します。また時間の入力にはコロン、 日付の入力にはダッシュも自動入力と なります。
最後に入力した文字を消去する	
入力した文字を全て消去する	Clear
入力画面を終了し、 変更箇所を保存する	Done

### 1.3 メニューの概要

この項では、主要な画面の概要について解説しています。

Main Menu 画面 システムにログインすると、Main Menu 画面が現れます。この画面は Veriti<sup>TM</sup> サーマルサイクラーを操作する初期画面になります。日付、時間、現在のユー ザー名が画面の下段に表示されます。

Main Menu						
Browse / New Methods	Settings Menu	Tools Menu	My Account			
Shortcut 1	Shortcut 2	Shortcut 3	Shortcut 4			
Shortcut 5	Shortcut 6	Shortcut 7	Shortcut 8			
Log Off 2007-01-1	16   3:42 PM	User: greg	or ?			

Veriti™ サーマルサイクラーの Main Menu 画面

Main Menu の各ボタンは:

- Browse/New Methods:このボタンから使用するすべてのメソッドファ イルを閲覧・作成することができます。
- Settings Menu: このボタンから装置の様々な初期設定を行なえます。アドミニストレーター権限でアクセスする場合は、この画面から Admin Menu 画面を表示することができます。
- Tools Menu: このボタンから装置の診断テストや Tm 計算機能などのツ ールを表示できます。
- My Account: このボタンからアカウント情報の閲覧・編集ができます。
   装置がログインを必要としていない設定の場合はこのボタンは表示されません。
- Shortcut:このボタンは、よく使う機能を Main Menu の画面上に登録 することができます。
- Log Off: このボタンで装置からログオフすることができます。但し、 ラン実行中は終了するまでログオフできません。
- (?):このボタンからヘルプ画面を表示します。

Settings Menu この画面から装置の様々な初期設定を行います。装置にログインした"User" 画面 または"Administrator"のみがこのメニューにアクセスできます。

Settings Menu						
Set Idle Temperatures	Set Time-Outs	Set Default Method	Configure Printer			
Configure Shortcuts	About the Instrument	Manage Folders	Admin Menu			
			To Previous Menu			
	2007-03-10	11:24 AM	?			

Settings Menu 画面。Admin Menu ボタンが表示されている。

### Admin Menu この画面は、ユーザーアカウントの操作、装置のセキュリティ設定、ランメソ 画面 ッドやセッティングのバックアップ、ファームウェアのアップグレード等の機 能を設定します。ログインが必要な設定の場合は、"Administrator"のみがこ のメニューにアクセスできます。



Admin Menu 画面。Administrator 権限があるユーザーのみ利用可能。

Tools Menu 画面 この画面から様々なツールや装置のテストにアクセスすることができます。



Tools Menu 画面。この画面はすべてのユーザーで利用可能。

Tools Menu の機能に関する詳細な情報は、英文 User Guide をご参照下さい。

## 1.4 ランメソッドについて

**ランメソッド** ランメソッドは、"どのような温度条件で装置を運転するか"を定義したもので、 とは 装置のファームウェアに保存されます。

**デフォルトで** Veriti<sup>™</sup> サーマルサイクラーのファーウェアには予め 10 個のランメソッドが 保存されている 保存されています。 ランメソッド

- AmpliCycle<sup>®</sup> Seq Run Method
- AmpliTaq Gold<sup>®</sup> DNA Polymerase Run Method
- BigDye<sup>®</sup> Standard Run Method
- BigDye<sup>®</sup> Fast Run Method
- Fast PCR Run Method
- General PCR Run Method
- LMS2 Run Method
- Time Release Run Method
- Touchdown PCR Run Method
- XL PCR Run Method

各ランメソッドは Shared フォルダに保存されています。これらランメソッド は異なるランメソッドとして編集・コピー・ランを行なうことができます。デ フォルトのランメソッドを消去することはできません。

### ランメソッドの 《ランメソッドの閲覧方法》

閲覧

 ランメソッド選択画面より、閲覧するメソッドをタッチして選択し、次 に View/Edit ボタンをタッチします。 ランメソッドが表示されます。



### !重要! "Limited User"または"Guest"としてログインしている場合 は、ランメソッドの変更、ラン実行は可能ですが、変更したファイルを保 存することはできません。

- 2. 閲覧する:
  - ランメソッド中の表示されていないステップを見るには、ランメソッドの端に表示されている矢印ボタンをタッチします。
  - 各ステップのランプレートを見るには、Options をタッチした後に Ramp Rate を選択します。画面上では各ステップの時間のフィールド に対応するランプレートが表示されます。
- 3. 閲覧が終了したら、×ボタンをタッチしてランメソッド選択画面に戻り ます。



### **ランメソッドの**現在のフォルダ内のランメソッドを見るには

管理

Main Menu 画面から Browse/New Methods をタッチします。
 ランメソッド選択画面が現れ、上段のタイトルには含まれるランメソッドの数がカッコ内に表示されます。

Run Method			Folder		Last	Used 🔻	
LMS2*			Shared		2007-	02-15	
Touchdown_P	CR*		Shared		2007-	02-15	
AmpliTaq_Goli	d*		Shared		2007-	02-14	Page 1/3
Time_Release			Shared		2007-	02-14	
XL_PCR*			Shared		2007-	02-14	<b>V</b>
		a de la		<u>ال</u>		<b>m</b>	Selected
Start Run	New	View/Edit	Сору	Prin	t	Delete	,
Touc to p	h a run met erform an a	hod to select i action. Touch a	t, then touch a column title	any of to sort	the bu the ta	ttons ble.	(?

ランメソッド選択画面。最近アクセスしたランメソッドが表示されます。

《ランメソッドの選択》

- 1. ランメソッドがすでにランメソッド選択画面で選択されている場合:
  - a. 選択されている行をタッチして、選択を解除します。
  - b. 目的のランメソッドの行をタッチして、選択します。
- 2. 1 つ以上のランメソッドを選択する場合は、各メソッドをタッチして選 択してください。

ランメソッドのリスト表示で操作可能な機能について下表に示します。

ランメソッドの管理

操作	使用するキー
ランメソッドを異なる基準で並べ替	ランメソッドの並べ替えをしたいカラムタイ
える	トルをタッチします。
	もう一度タイトルをタッチすると、逆の順番に
	並べ替えられます。
フォルダ内にはあるが、画面には表	<b>、</b> をタッチします。
示されていないランメソッドを見る	
異なるフォルダ内のランメソッドを	
見る	ストが表示されるので、閲覧したいフォルダを
	タッチして選択します。
ランを開始する	ランメソッドのリストで、ランメソッドをタッ
	チして選択した後、Start Run をタッチします。
	(p38 の"ランの実行"を参照してください)

操作	使用するキー
新しいランメソッドを作成する	New をタッチします。(p29 "新しいランメソ
	ッドの作成"を参照してください)
ランメソッドの閲覧または編集を	ランメソッドのリストから目的のランメソッ
する	ドをタッチして選択した後、View/Edit をタッ
	チします。(p31"ランメソッドの編集"を参照
	してください)
ランメソッドをコピーする	ランメソッドのリストから目的のランメソッ
	ドをタッチして選択した後、Copy をタッチし
	ます。(p14 "ランメソッドのコピー"を参照し
	てください)
ランメソッドを印刷する	ランメソッドのリストから目的のランメソッ
	ドをタッチして選択した後、 <b>Print</b> をタッチし
	ます。
ランメソッドを消去する	ランメソッドのリストから目的のランメソッ
	ドをタッチして選択した後、 <b>Delete</b> をタッチし
	ます。(p15 "ランメソッドの消去"を参照して
	ください)
ランメソッドをデフォルトとして	p15"ランメソッドをデフォルトとして設定"
設定する	を参照してください。
ランメソッドをバックアップする	p16 "ランメソッドのバックアップとその他の
	設定"を参照してください。
以前バックアップしたランメソッ	p17 "ランメソッドやその他の設定の回復・転
ドを再度保存する	送"を参照してください。

ランメソッドの管理(つづき)

ランメソッドの

コピー

《ランメソッドのコピー方法》

- 1. "User"または"Administrator"でログインをします。
- 2. ランメソッド選択画面で、コピーをするランメソッドを確認します。
- 3. 目的のランメソッドをタッチした後、**Copy**をタッチします。
- 4. Save Method 画面において
  - a. コピーするランメソッドが1つの場合は、Name をタッチしてメ ソッドの新しい名前をつけます。そのまま同じフォルダに保存す るか、Folder をタッチしてから異なるフォルダに保存します。 コピーするランメソッドが複数の場合は、Folder をタッチして ランメソッドを異なるフォルダに保存します。
  - b. 必要なら、**Reaction Volume**, **Cover Temperature**, **Notes** などをタ ッチして選択し、数値の変更を行ないます。また、このメソッド を実行するときにこれらの値を変更することもできます。
  - c. Save & Edit をタッチした後、OK を選択します。

NOTE ログインが必要な設定になっている場合は、"User"または"Administrator"レベルの権 限がないとコピー操作はできません。ログイン設定やセキュリティに関する内容については、p47 "セキュリティおよびユーザーアカウントの設定"をご参考下さい。

### **ランメソッドの** 《ランメソッドの消去方法》

消去

- 1. "User"または"Administrator"でログインをします。
- 2. ランメソッド選択画面で、消去するランメソッドを確認します。
- 3. 目的のランメソッドをタッチした後、Delete をタッチします。
- 4. ダイアログボックスが表示され、消去してよいかの確認をします。
- 5. メッセージボックスの Yes をタッチするとランメソッドは消去されます。

NOTE ログインが必要な設定になっている場合は、"User"または"Administrator"レベルの権 限がないとランメソッドの消去はできません。ログイン設定やセキュリティに関する内容について は、p47 "セキュリティおよびユーザーアカウントの設定"をご参考下さい。

ランメソッドを
 デフォルト
 として設定
 フォルト
 として設定
 フォルト設定をしていないと
 Vyドを使用します。

### !重要! ログインが必要な設定になっている場合、ランメソッドをデフ オルトとして設定するには、"User"または"Administrator"でログイン する必要があります。Shared Folderに保存されているランメソッドのみが、 デフォルトのランメソッドとして設定できます。

《デフォルトのランメソッドの設定方法》

- 1. "User" または "Administrator" でログインをします。
- 2. Main Menu 画面より、Settings Menu をタッチします。
- Settings Menu 画面で、Set Default Method をタッチします。
   デフォルトランメソッド設定画面には、Shared Folder にあるランメソッドのみが表示されます。
- デフォルトランメソッド設定画面より、目的のランメソッドをタッチした後、Make Default をタッチします。メッセージボックスが表示されるので内容を確認した後、OK をタッチします。 Settings Menu 画面が表示され、選択したランメソッドがデフォルトとして設定されます。

**ランメソッド** 装置のランメソッドとその他の設定は、USB ドライブにバックアップすること **及び各設定の** ができます。すべてのランメソッド、フォルダ、ユーザー情報、その他装置の **バックアップ** 設定などが保存されます。

> NOTE: ログインが必要な設定になっている場合は、"Administrator"レベルの権限がないとラ ンメソッドやセッティングのバックアップはできません。ログイン設定やセキュリティに関する内 容については、p47 "セキュリティおよびユーザーアカウントの設定"をご参考下さい。

《ランメソッドやその他の設定の回復・転送方法》

1. USB ドライブを装置正面の USB ポートに挿入します。



- 2. "Administrator"として装置にログインをします
- 3. Main Menu 画面から Settings Menu、Admin Menu の順番にタッチしま す。
- 4. Admin Menu 画面から Back Up Methods & Settings をタッチします。 Back Up Methods & Settings 画面が開きます。

Back Up Methods and Settings		
Back Up		
Insert a USB drive, then touch Back Up to save all run methods and instrument settings to the memory stick.	?	

5. Back Up をタッチします。

バックアップファイルがすでに存在する場合は、上書きを確認する画面 が表示されます。 装置は USB へのすべてのランメソッド、フォルダ、ユーザー情報やそ の他の情報のコピーが完了すると、バックアップの完了を示すメッセー

の他の情報のコピーが完了すると、バックアップの完了を示すメッセジボックスが表示されます。

- 6. OK をタッチし、× をタッチすると Admin Menu 画面に戻ります。
- 7. USB ドライブを取り外し、安全な場所に保管します。

ランメソッドや
 USBドライブに保存されているランメソッドやその他の設定を、装置に回復さ
 その他の設定の
 せることができます。また、このデータ転送の機能は、データ元以外の装置に
 回復・転送
 もデータを転送・保存することができます。

NOTE: ログインが必要な設定になっている場合は、"User"または"Administrator"レベルの 権限がないと、保存してあるランメソッドやその他の設定を装置に転送することはできません。ロ グイン設定やセキュリティに関する内容については、p47"セキュリティおよびユーザーアカウン トの設定"をご参考下さい。

《ランメソッドやその他の設定の回復・転送方法》

- 1. ランメソッドやその他の設定を保存した USB ドライブを装置正面の USB ポートに挿入します。
- 2. "User"または"Administrator"として装置にログインをします。
- 3. Main Menu 画面から Settings Menu、About the Instrument の順番にタッ チします。



- About the Veriti<sup>™</sup> Thermal Cycler 画面で、Restore Method & Settings をタッチするとメッセージボックスが現れるので、Yes をタッチします。 装置はすべてのランメソッド、フォルダ、ユーザー情報やその他の情報 を USB ドライブから装置にコピーを開始し、完了するとメッセージボ ックスが現れます。
- 5. USB ドライブを取り外し、安全な場所に保管します。
- 6. 装置背面のスイッチを使用して装置の電源を切ります。
- 7. 装置の電源を入れて、再起動をさせます。

!重要! データ元以外の装置にバックアップデータを転送した場合、その装置がネットワーク上に存在しているのであれば、装置の名称を変更する必要があります(英文ユーザーガイドの P2-8 Setting the instrument name"を参照してください)。

## 1.5 ランフォルダについて

ランメソッドは、いずれかのフォルダに保存されます。そして出荷時のデフォ ルト状態では3つのフォルダが選択できます。

- Default-特にフォルダの指定をしない限り、ランメソッドはこのフォ ルダに保存されます。ランメソッドはファイルを作成したユーザーで あれば、閲覧、編集、実行、消去をすることができます。デフォルト フォルダは青いボタンです。
- Shared-すべてのユーザーが、ランメソッドのアクセス、編集、実行、 消去を行なえます。フォルダ選択画面において Shared フォルダは青い ボタンで表示されます。
- USB-このフォルダはUSBドライブに保存されたすべてのランメソッドを表示します。USBドライブが装置に接続されている場合は青いボタンで、接続されていない場合は、グレーで表示されます。

カスタムフォルダは9個まで作成することができます。カスタム設定したフォ ルダでは、ランメソッドを作成したユーザーであれば、閲覧、編集、実行、消 去をすることができます。フォルダ選択画面で、カスタムフォルダは黄色で表 示されます。 ランメソッドを保存するときに、新しいフォルダを作成することができます (p30の操作ステップ6を参照)。

- **カスタムフォルダ** カスタムフォルダは消去することができます。カスタムフォルダを消去すると、 の消去 フォルダに含まれるすべてのランメソッドも消去されます。
  - 1. Main Menu 画面から Settings Menu をタッチします。
  - 2. Settings Menu 画面の Manage Folders をタッチします。
  - 3. Manage Folders 画面で、消去したいフォルダのチェックボックスにタ ッチして選択した後、**Delete** をタッチします。
  - 4. ダイアログボックスが現れますので、内容を確認して Yes、その後 OK をタッチします。
  - 5. 引続き別のフォルダを消去する場合は、上記ステップ3と4を繰り返し ます。
  - 6. フォルダの消去が完了したら、 × をタッチして Settings Menu 画面に 戻ります。

### カスタムフォルダ

カスタムフォルダの名称は変更することができます。

の名称変更

- 1. Main Menu 画面から Settings Menu をタッチします。
- 2. Settings Menu 画面の Manage Folders をタッチします。
- 3. Manage Folders 画面で、名称変更をしたいフォルダのチェックボック スにタッチして選択した後、**Rename** をタッチします。
- 4. フォルダの新しい名称を最大 12 文字の範囲で入力をした後、**Done** をタ ッチすると Manage Folders 画面に戻ります。
- 5. 引続き別のフォルダの名称を変更する場合は、上記ステップ3と4を繰り返します。
- 6. フォルダの消去が完了したら、 × をタッチして Settings Menu 画面に 戻ります。

## 2. ランの実行

このセクションでは下記について説明しています。

- 2.1 消耗品の選択
- 2.2 サンプルを装置にセットする
- 2.3 新しいランメソッドの作成
- 2.4 詳細なメソッドの編集
- 2.5 ランの実行
- 2.6 ランのモニタリング
- 2.7 ランレポートの閲覧と印刷
- 2.8 装置からサンプルを取り出す
- 2.9 サイクルパフォーマンステストの実行

### 2.1 消耗品の選択

このセクションでは、Veriti<sup>TM</sup>サーマルサイクラーで使用する最適なチューブ、 プレート、サンプルトレーについて解説しています。

!重要! サンプルチューブを使用して Veriti<sup>™</sup> サーマルサイクラーを 運転する場合は、必ずトレーまたはトレー/リテーナーを使用してください

**チューブの選択** 次の3種類のサンプルチューブが利用できます。 以下のVeriti<sup>™</sup> サーマルサイクラーに使用する3種類のチューブの組み合わせ を参照して、消耗品のご用意をお願いします。

キャップ付シングルチューブ:

MicroAmp® Fast Reaction Tube with Cap(0.1mL) または MicroAmp® Reaction Tube with Cap(0.2mL)



MicroAmp<sup>®</sup> Reaction Tube  $(0.2mL) \succeq$ MicroAmp<sup>TM</sup> 8-Cap strip



8連チューブと8連キャップ:

MicroAmp<sup>™</sup> Fast 8-Cap strip (0.1mL)と MicroAmp<sup>™</sup> 8-Cap strip または MicroAmp<sup>™</sup> 8-Tube Strip(0.2mL) と MicroAmp<sup>™</sup> 8-Cap strip



### サンプルトレー とプレートの 組み合わせ

### **サンプルトレー** サンプルトレーとプレート組み合わせを以下の表に示します。

! 重要! サンプルをチューブまたはプレートに分注する際は MicroAmp<sup>™</sup> Splash Free 96-Well Base をご使用ください。但し、この Base はサンプルブロックにはセットしません。

Run Vessel	Compatible Component	Illustration
MicroAmp <sup>™</sup> Fast Optical 96-Well Reaction Plate (0.1-mL) MicroAmp <sup>™</sup> Optical 96-Well Reaction Plate (0.2-mL)	MicroAmp <sup>™</sup> Caps, 8-Cap Strip or MicroAmp <sup>™</sup> Clear Adhesive Film (not shown)	MicroAmp <sup>™</sup> 8-Cap Strip MicroAmp <sup>™</sup> Fast Optical 96-Well Reaction Plate (0.1-mL) or MicroAmp <sup>™</sup> Optical 96-Well Reaction Plate (0.2-mL) MicroAmp <sup>™</sup> Splash Free 96-Well Base
MicroAmp <sup>®</sup> Fast Reaction Tube with Cap (0.1-mL) MicroAmp <sup>®</sup> Reaction Tube with Cap (0.2-mL)	96-Well Tray for VeriFlex™ Blocks	MicroAmp <sup>®</sup> Fast Reaction Tube with Cap (0.1-mL) or MicroAmp <sup>®</sup> Reaction Tube with Cap (0.2-mL) MicroAmp <sup>™</sup> 96-Well Tray for VeriFlex <sup>™</sup> Blocks MicroAmp <sup>™</sup> Splash Free 96-Well Base
MicroAmp <sup>™</sup> 8-Tube Strip (0.2-mL) <i>or</i> MicroAmp <sup>®</sup> Reaction Tube without Cap (0.2-mL)	MicroAmp <sup>™</sup> 8-Cap Strip and MicroAmp <sup>™</sup> 96-Well Tray/Retainer Set for Veriti <sup>™</sup> Systems	MicroAmp <sup>™</sup> 8-Cap Strip MicroAmp <sup>™</sup> 96-Well Retainer MicroAmp <sup>™</sup> 8-Tube Strip (0.2-mL) or MicroAmp <sup>®</sup> Reaction Tube without Cap (0.2-mL) MicroAmp <sup>™</sup> 96-Well Tray for Veriti <sup>™</sup> Systems MicroAmp <sup>™</sup> Splash Free 96-Well Base

Part Description	96-Well (0.1-mL)	96-Well (0.2-mL)	Part Number		
96-Well Plates					
MicroAmp <sup>™</sup> Fast 96-Well Reaction Plate (0.1 mL)	x		4346907		
MicroAmp <sup>™</sup> Fast Optical 96-Well Reaction Plate	x		4346906		
	x		4366932		
MicroAmp™ Optical 96-Well Reaction Plate		х	N8010560		
		х	4316813		
MicroAmp <sup>™</sup> Optical 96-Well Reaction Plate with		х	4306737		
Barcode		х	4326659		
MicroAmp™ Optical 96-Well Reaction Plate with Barcode and Optical Caps		х	403012		
MicroAmp™ Optical 96-Well Reaction Plate with Barcode and Optical Adhesive Films		х	4314320		
12-Well Strips					
MicroAmp™ 12-Cap Strip	x	х	N8010534		
	x	х	N8011534		
MicroAmp™ 12-Cap Strip, Assorted Colors	x	х	N8010834		
8-Well Strips					
MicroAmp™ Fast 8-Tube Strip (0.1 mL)	х		4358293		
MicroAmp™ 8-Tube Strip (0.2 mL)		х	N8010580		
MicroAmp™ 8-Tube Strip, Assorted Colors (0.2 mL)		х	N8010838		
MicroAmp™ Optical 8-Tube Strip (0.2 mL)		х	4316567		
MicroAmp™ 8-Cap Strip	x	х	N8010535		
	x	х	N8011535		
MicroAmp™ 8-Cap Strip, Assorted Colors	x	х	N8010835		
MicroAmp™ Optical 8-Cap Strip	x	х	4323032		
Single Tubes	•				
MicroAmp® Fast Reaction Tube with Cap (0.1 mL)	х		4358297		

Veriti<sup>TM</sup> サーマルサイクラー用 消耗品

Part Description	96-Well (0.1-mL)	96-Well (0.2-mL)	Part Number
MicroAmp® Reaction Tube with Cap (0.2 mL)		x	N8010540
		х	N8011540
MicroAmp® Reaction Tube with Cap, Assorted Colors (0.2 mL)		x	N8010840
MicroAmp® Reaction Tube with Cap, Autoclaved (0.2 mL)		x	N8010612
MicroAmp® Reaction Tube without Cap (0.2 mL)		x	N8010533
		х	N8011533
MicroAmp® Reaction Tube without Cap, Assorted Colors (0.2 mL)		x	N8010833
MicroAmp® Optical Tube without Cap (0.2 mL)		х	N8010933
Seals & Covers			
MicroAmp™ Clear Adhesive Film	x	x	4306311
MicroAmp™ Optical Adhesive Film	x	x	4360954
	x	x	4311971
MicroAmp <sup>™</sup> Optical Adhesive Film Kit	x	х	4313663
Reaction Trays			
MicroAmp™ 96-Well Tray for VeriFlex™ Blocks	x	x	4379983
MicroAmp <sup>™</sup> 96-Well Tray/Retainer Set for Veriti <sup>™</sup> Systems		x	4381850

Veriti™ サーマルサイクラー用 消耗品 (続き)

その他のサンプル調製用備品

次の消耗品はサンプル調製時や Veriti<sup>TM</sup> サーマルサイクラー稼動時に使用する と便利です。

Product	Part Number
MicroAmp™ 96-Well Base	N8010531
MicroAmp™ Splash Free 96-Well Base	4312063
MicroAmp™ Cap Installing Tool (Handle)	4330015
MicroAmp™ Multi-Removal Tool	4313950
MicroAmp <sup>™</sup> Adhesive Film Applicator	4333183
MicroAmp™ Centrifuge Adapter	N8013822

## 2.2 サンプルを装置にセットする

このセクションでは以下のサンプルチューブを装置にセットする方法について 解説しています。

- キャップ付きチューブ
- 96-well プレートタイプのチューブ
- キャップなしチューブ(8連チューブ、キャップなしシングルチューブ)

サンプルブロックウェルやチューブ内に、蒸発防止としてミネラルオイルやグ リセリンを使用する必要はありません。MicroAmp® Reaction Tube は、ブロッ クウェルに密着し、装置のヒーターカバーはどの位置のチューブに対しても均 等に加温・加圧するため、チューブ内での結露を排除するように設計・製造さ れています。

### Veriti サーマルサイクラー: 0.2mL ブロックの場合

キャップ付き チューブの	1.	96 ウェルトレーを Splash Free 96-Well Base の上にセットします。
セット	2.	トレーにキャップ付きチューブをセットします。
	3.	サンプルをチューブに分注します。
	4.	チューブにキャップをします。
	5.	p28 の"サンプルトレーまたはプレートをサンプルブロックにセット" を参照し、装置にセットをします。
96-well プレート のセット	1.	96-well プレートを Splash Free 96-Well Base の上にセットします。
	2.	サンプルをチューブに分注します。
	3.	次のいずれかを使用してふたをします。
		• MicroAmp <sup>TM</sup> 8-Cap strip
		• MicroAmp <sup>TM</sup> 12-Cap strip
		• MicroAmp <sup>TM</sup> Optical 8-Cap strip
		• MicroAmp <sup>TM</sup> Clear Adhesive Film

- MicroAmp<sup>TM</sup> Optical Adhesive Film
- p28の"サンプルトレーまたはプレートをサンプルブロックにセット" を参照し、装置にセットをします。

#### キャップなし チューブの セット 1. 96 ウェルトレー/リテーナーのトレーを Splash Free 96-Well Base の上 にセットします。

- 2. シングルまたは8連チューブをトレーにセットします。
  - 3. リテーナーをチューブの上からセットします。
  - 4. サンプルをチューブに分注します。
- 5. MicroAmp<sup>™</sup> 8-Cap Strip でチューブにキャップをします。
- p28の"サンプルトレーまたはプレートをサンプルブロックにセット" を参照し、装置にセットをします。

### Veriti Fast サーマルサイクラー: 0.1mL ブロックの場合

キャップ付き

- チューブの
- 1. 96 ウェルトレーを Splash Free 96 Well Base の上にセットします。
  - **セット** 2. トレーに Fast キャップ付きチューブをセットします。
    - 3. サンプルをチューブに分注します。
    - 4. チューブにキャップをします。
    - p28の"サンプルトレーまたはプレートをサンプルブロックにセット" を参照し、装置にセットをします。

96-well Fast プレートを Splash Free 96-Well Base の上にセットします。

96-well Fast

プレートの

セット

- 2. サンプルをチューブに分注します。
  - 3. 次のいずれかを使用してふたをします。
    - MicroAmp<sup>TM</sup> 8-Cap strip
    - MicroAmp<sup>TM</sup> 12-Cap strip
    - MicroAmp<sup>TM</sup> Optical 8-Cap strip
    - MicroAmp<sup>TM</sup> Clear Adhesive Film
    - MicroAmp<sup>TM</sup> Optical Adhesive Film
  - p28の"サンプルトレーまたはプレートをサンプルブロックにセット" を参照し、装置にセットをします。

ast

Fast 8 連チューブ のセット

- 1. 96 ウェルトレーを Splash Free 96-Well Base の上にセットします。
  - 2. トレーに Fast 8 連チューブをセットします。
  - 3. サンプルをチューブに分注します。
  - 4. MicroAmp<sup>™</sup> 8-Cap Strip でチューブにキャップをします。
  - 5. p28 の "サンプルトレーまたはプレートをサンプルブロックにセット" を参照し、装置にセットをします。

**サンプルトレー** サンプルブロックにサンプルトレー またはプレートを せ〈アセンブリ〉においても同様 サンプルブロック

にセット

- サンプルブロックにサンプルトレーをセットする手順は、以下のいずれの組合 せ〈アセンブリ〉においても同様です。
  - トレー/リテーナーを使用したチューブアセンブリ
  - トレーのみでチューブを保持するアセンブリ
  - 96 ウェルプレートによるアセンブリ
  - 1. サンプルチューブアセンブリまたは 96-well プレートを Splash Free 96-Well Base から取り外し、装置のサンプルブロックにセットします。
  - 2. セットするアセンブリまたは 96-well プレートの向きは、以下に示すようにウェルポジション A1 が左上になるようにします。





!重要! ベースはサンプルブロックにセットしません。

3. 装置のヒーターカバーを閉めます。

## 2.3 新しいランメソッドの作成

新規作成

**ランメソッドの** 《ランメソッドの新規作成方法》

- 1. Main Menu 画面から Browse/New Methods をタッチします。
  - ランメソッド選択画面から New をタッチします。 デフォルトのランメソッドが Edit Run Method 画面に表示されます(デ フォルトのランメソッド変更したい場合には、p 15 の "ランメソッドを デフォルトとして設定"を参照)

Edit P	un Methoo	I: LMS2		×
Stage 1 × 1	Stage 2 × 10		Stage 3 × 20	
95.0	94.0	55.0         72.0           0:15         0:30	89.0 0:15	55.0
Step 1	Step 1	Step 2 Step 3	Step 1	Step 2
	d <sup>i</sup> o			0+
Run Touch a or temp or Ve	Add stage or a ste erature to edi riFlex steps, t	Delete p to insert a stage or st t it. Touch Options to c o show ramp rates or to	Options zep. Touch a time reate AutoDelta add a pause.	Save

- 3. デフォルト表示されたランメソッドから目的の設定条件に編集します。
  - ステージの追加 (p31 の "ステージの追加"を参照)
  - ステップの追加 (p32 の "ステップの追加"を参照)
  - VeriFlex ブロックに異なる温度プログラムのステップを追加 (p34 の "VeriFlex ブロックで異なる温度プログラムを設定"を参照)
  - AutoDelta ステップの追加 (p35 の "AutoDelta ステップの設定"を 参照)
  - ポーズ設定の追加 (p37 の "ポーズの設定"を参照)
  - ステップまたはステージの消去(p37の"ステップまたはステージの 消去"を参照)

4. Save & Exit をタッチすると、Save Run Method 画面が表示されます。

Save Run Methoo	
Run Method:	LMS2
Folder:	Shared
Reaction Volume: Notes:	10     Cover Temperature:     105.0       This method used for gene expression al <sup>a</sup> ssay.
	Save & Exit
Enter a name, rea for this run me When you are	ction volume, and choose a folder thod. Notes can also be added. finished, touch 'Save & Exit'.

- 5. Run Method のフィールドをタッチして、ランメソッドの名称を入力しま す。名称はスペースを含まない最大 48 文字まで使用することができま す。
- 6. Folder のフィールドをタッチしてランメソッドを保存するフォルダを選 択します。
  - すべてのユーザーがアクセスできるように保存するには、Sharedを選択してください。
  - 新しいフォルダを作成するには、リスト下段の New Folder を選択し、 スペースを含まない最大 12 文字の名称を入力した後、Done をタッチ してください。
  - ランメソッドを USB ドライブに保存するには、USB をタッチしてく ださい(USB ドライブが装置に装着されているときのみ利用可能で す)。
- 7. Reaction Volume をタッチし、反応液量を入力します(このランメソッド を運転するときに、改めて反応液量を入力できます) v
- 8. Cover Temperature をタッチし、装置のヒーターカバーの温度を設定しま す(このランメソッドを運転するときに、改めて設定温度を入力できま す)。
- 9. Notes をタッチすると、このランメソッドに関するメモを入力できます。
- 10. 入力が完了したら Save & Exit をタッチします。 ランメソッド選択画面に戻ります。

## 2.4 ランメソッドの編集

このセクションでは以下に記すランメソッドの編集方法について記します

- ステージの追加 (p31)
- ステップの追加 (p32)
- 温度と時間の変更(p33)
- 温度の移行速度(ランプレート)の変更(p33)
- VeriFlex ブロックで異なる温度プログラムを設定(p34)
- AutoDelta ステップの設定 (p35)
- ポーズの追加 (p37)
- ステージまたはステップの消去 (p37)
- **ステージの追加** ランメソッドのステージは、2つのタイプに大別されます。
  - サイクルステージ: サイクルステージは、設定温度を指定されたサイクル数繰り返します。一般的なサイクルステージのステップは、テンプレート DNA の変性、プライマーのアニーリング、プライマーの伸長の3ステップからなります。
  - ホールドステージ: ホールドステージは、1サイクルのみの反応です。

《ステージの追加方法》

1. Edit Run Method 画面でステージ上部のステージ名をタッチします。ス テージが選択されるとステージ名の近傍エリアが赤になります。

Edit F	Run Method: I	_MS2		×
Stage 1 × 1 95.0	Stage 2           ×         10           94.0	55.0 72.0	Stage 3 × 20 89.0	55.0
12:00	0:15	0:15 0:30 Step 2 Step 3	0:15	0:15 Step 2
	o <sup>t</sup> o		30 ▲	0+
Run Touch a or temp or Ve	Add stage or a step t perature to edit it. riFlex steps, to s	Delete o insert a stage or sta . Touch Options to cr how ramp rates or to	Options ep. Touch a time eate AutoDelta add a pause.	Save ?

 Add をタッチします。 選択したステージのコピーがランメソッドに追加されます。

- 3. サイクルステージには、必要なサイクル数を設定します。
  - a. ステージをタッチで選択します。
  - b. サイクル数の (ステージ名の下にある) 入力フィールドをタッチ します。

	Edit Ru	n Method	: LMS2				
Sta ×	age 1 1	Stage 2 × 10	<		Stage 3 × 20		
<b>9</b>   <b>1</b>	15.0 [ 2:00 [	94.0 0:15	55.0 0:15	72.0	89.0 0:15	<u>55.0</u> 0:15	
St	ep 1	Step 1	Step 2	Step 3	Step 1	Step 2	
(	1	2	з	4	5		
(	6	7	8	9	0	Do	ne

c. サイクル数を入力した後、**Done**をタッチします。

4. 温度や時間を変更する場合は、p33の"温度と時間の変更"を参照して ください。

### **ステップの追加** 《ステップの追加方法》

1. Edit Run Method 画面で、選択するステップのカラム内をタッチしま す。ステップが選択されると、カラムは赤くなります。



 Add をタッチします。 選択したステップのコピーがランメソッドに追加されます。

Edit R	un Methoo	I: LMS2		×
Stage 1	Stage 2			Stage 3
× 1 95.0	94.0 0:15	94.0 0:15 0:15	72.0	89.0 0:15
Step 1	Step 1	Step 2 Step 3	Step 4	Step 1
	o <sup>t</sup> o			0+
Run Touch a or temp or Ve	Add stage or a ste erature to edi riFlex steps, t	Delete p to insert a stage or sta t it. Touch Options to cr o show ramp rates or to	Options ep. Touch a time eate AutoDelta add a pause.	Save ?

- 3. 各ステップの温度や時間を変更する場合は、p33の"温度と時間の変更" を参照してください
- 温度と時間の変更
- 1. 変更したい温度または時間のフィールドをタッチし、温度・時間の数値 を入力します。
- 2. その他の温度・時間パラメーターを変更したい場合は、編集画面を閉 じずにそのまま変更・設定を行ないます。
  - a. 変更するフィールドをタッチします。
  - b. 編集画面で数値をタッチ入力します。
  - c. その他のフィールドを変更する場合は、上記操作を繰返します。
- 3. Done をタッチすると編集画面が閉じます。

**温度の移行速度の** ステップ間の温度の移行速度(ランプレート)を各ステップごとに設定するこ 変更 とができます。デフォルトのランプレートは、すべて 100%です。

- 1. 変更したいランプの右側のステップをタッチします。
- 2. Options をタッチし、Ramp Rate を選びます。画面上の時間のパラメー タがランプレートの%表示に変わります。
- 3. ランプレートの入力フィールドをタッチし、任意のランプレートの数値 を入力します。ランプレートの入力範囲は 0.1~100%です。
- Done をタッチすると編集画面が閉じます。
   ランプレートを 100%以外に設定すると変更・設定されたステップの隣 に配のシンボルが表示されます。

VeriFlex ブロックで 異なる温度プログラム を設定 Veriti<sup>TM</sup> サーマルサイクラーのサンプルブロックは、6つの温度ゾーンをもち、 各ゾーンは独立の温度設定ができます。またすべてのブロックを単一のゾーン として運転することもできます。デフォルトでは単一温度ゾーンとして設定を します。VeriFlex<sup>TM</sup>ブロック上で異なる温度設定で運転するステップを、ここ では"VeriFlex<sup>TM</sup>ステップ"と呼びます。この機能は、新たな PCR 条件の最 適化にたいへん有用です。

《VeriFlex<sup>TM</sup>ブロックで異なる温度のステップを作成》

1. 温度設定をしたいステップをタッチして選択した後、**Options** をタッチ するとオプションメニューが開きます。

Edit P	Run Methoo	I: LMS2		×		
Stage 1 × 1	Stage 2 × 10		Stage 3 × 20			
95.0	94.0	55.0         72.0           0:15         0:30	89.0 0:15	55.0 0:15		
Step 1	Step 1	Step 2 Step 3	Step 1	Step 2		
	o <sup>t</sup> o			<b>O</b> •		
Run	Add	Delete	Options	Save		
Touch a stage or a step to insert a stage or step. Touch a time or temperature to edit it. Touch Options to create AutoDelta or VeriFlex steps, to show ramp rates or to add a pause.						

2. オプションメニュから VeriFlex Step をタッチすると、VeriFlex ステップ作成画面が開きます。

Edit Ru	n Method	: LMS2			
	C	reate a Ve	eriFlex St	ер	?
55.0	56.0	57.0	58.0	59.0	60.0
1 – 2	3 - 4	5 - 6	7 - 8	9 - 10	11 - 12
*Tempera	ture differer	ice between	adjacent zo	ones <= 5.0	Apply Selected to All
1	2	3	4	5	
6	7	8	9	0	Cancel Done

 温度フィールドをタッチし、目的の温度を入力します。複数のゾーンを 同じ温度に設定することもできます。ただ異なる温度を設定する場合、 隣接するゾーン間の温度差は0.1℃から5.0℃の範囲で設定できます。 また最高設定温度は99.9℃です。 6つの各々のゾーンについて、上記ステップ3を繰り返してください。
 入力が完了したら Done をタッチします。

*NOTE: すべてのフィールドに対して、同一温度を入力する場合は、まずひとつのフィールドに 目的の温度を入力した後、Apply Selected to All をタッチします。* 

### AutoDelta ステップ AutoDelta ステップ機能を使用すると、サイクルステージにおける温度や時間 の設定 のパラメータをサイクル毎に自動増減することができます。 AutoDelta ステップの応用例として、タッチダウン PCR があります。この PCR テクニックは初期のサイクルにおいて、アニーリング温度を高く設定し、サイ クルが進むにしたがいアニーリング温度をサイクル毎に徐々に低くしていく 技術です。このようなタッチダウン PCR を使用すると、PCR 産物の特異性を 高めることが可能になります。初期の高温条件ではプライマーとテンプレート の特異的アニーリングを助長させます。特異的に増幅したフラグメントは後半 に続く低温でのサイクルで有効にテンプレートとして PCR 増幅に利用される ため、低温におけるプライマーの非特異的な増幅を抑制します。

NOTE: AutoDelta ステップはサイクルステージでのみ設定が可能です。

《AutoDelta ステップの設定》

1. AutoDelta を設定したいステップをタッチした後、**Options** をタッチします。



2. AutoDelta を選択すると、AutoDelta ステップ設定画面が現れます。

Edit Run Method: LMS2	
Create an AutoDelta Step	?
Starting Cycle: 2	<u> </u>
Delta Temperature: + 0.0 deg C	
Delta Time: + 00:00 mm:ss	Remove AutoDelta Step
1 2 3 4 5	
6 7 8 9 0	Cancel Done

- 3. **Starting Cycle** をタッチして、AutoDelta ステップを開始するサイクル 番号を入力します。
- 4. 増減させるパラメーターを選択します。
  - 温度を増減したい場合: Delta Temperature をタッチして、サイクルごとに変更する温度を入力します。温度を低温側に変更する場合は、+(プラス)ボタンをタッチして-(マイナス)表示にします。+(プラス)表示の場合は、温度はサイクルごとに上昇します。
  - 時間を増減したい場合: Delta Time をタッチして、サイクルごとに変更する時間を入力します。時間をサイクルごとに減少させる場合は、+(プラス)ボタンをタッチして-(マイナス)表示にします。
     +(プラス)表示の場合、時間はサイクルごとに増加します。

NOTE: 温度の自動増減設定を行う場合、サイクル終了にともなう最終的な設定温度は 4.0℃以 上、99.9℃以下であることを確認してください。また減少させる時間の値は、従来設定してある保 持時間以下に設定していることを確認してください。

5. OK をタッチして設定を保存すると、ランメソッドを表示している画面 に戻ります。AutoDelta を設定したステップには AutoDelta のシンボ ル ▲が現れます。以下に示す図では、ステージ2のステップ2に AutoDelta の設定がされています。



**ポーズの追加** サイクルステージでは、任意のステップの前にポーズを設定することができます。

《ポーズの追加方法》

- 1. Edit Run Method 画面から、サイクルステージ内のポーズを設定した いステップをタッチして選択した後、**Option** をタッチします。
- 2. オプションメニューで Pause をタッチします。

Edit Ru	ın Method	: LMS2			
	Add	a Paus	e Before a	step	?
S	start first pau	se at cycl	le 1	of 10	
ı	<sup>2</sup> ause every	10	cycles for	00:10	mm:ss
					Remove Pause
1	2	3	4	5	
6	7	8	9	0	Cancel Done

- 3. 以下の項目を入力します。
  - ポーズを行なう最初のサイクル番号
  - ポーズの頻度
  - ポーズをする時間(分:秒)
- 4. Done をタッチしてポーズの設定を保存し、メソッド編集画面に戻りま す。ポーズ設定を行ったステップには ■というシンボルがメソッド上 に表示されます(以下の画面においてはステージ3のステップ3を参 照)。

Stage 2 × 10	Stage 3 × 20
<b>94.0</b>	60.0         72.0         89.0         55.0         72.0           0:15         0:30         0:15         0:30
Step 1	Step 2 Step 3 Step 1 Step 2 Step 3

**ステージまたは** 《ステージまたはステップの消去方法》

ステップの消去

- 1. Edit Run Method 画面で消去したいステージまたはステップをタッチ して選択します。
- 2. Delete、OK の順にタッチします。 ステージまたはステップはランメソッドから消去されます。

## 2.5 ランの実行

**ランを開始する前に** サンプル調製とサンプルのブロックへのセットに関しては、p26の"サンプル を装置にセットする"を参照してください。

### !重要! 2台以上の装置をネットワーク接続している場合、それら複数 台を同時に運転開始させることもできます。

《Main Menu 画面からラン開始》

- 1. 装置のカバーを閉めて、画面の Browse/New Methods をタッチします。
- 2. 使用するランメソッドをタッチして選択します。必要であれば、各パラ メータ設定が正しいか確認します。(p12 の "ランメソッドの閲覧"を 参照してください)。
- 3. Start Run をタッチすると、Run Parameters 画面が現れます。

Run Paramet	ers	×
Reaction Volume:	10	uL
Cover Temperature:	105.0	°C
Run ID:	200702140925	1
Notes:	This is a quick test for Joe's samples with a 5 ° minute vortex step.	
	Start Run Now Start	Multiple Runs
Touch to edit th	each field then use the keyboard e contents. When you are finished, touch Start Run Now.	?

- 4. Reaction Volume をタッチして反応液量を入力します。デフォルトの値 はランメソッドを作成・保存した時の値が表示されます 反応液量の範囲は装置のタイプによって異なります
  - Veriti<sup>™</sup> 96 ウェル Fast サーマルサイクラー: 5~40 µ L (10~30 µ L)\*
  - Veriti<sup>™</sup> 96 ウェルサーマルサイクラー: 10~100 µ L (10~80 µ L)\*

\* 上記の反応液量の範囲は、スクリーン上で入力可能な装置の温度管理能力範囲を示します。尚、 推奨する反応液量に関しては別途カッコ書きで併記しました。

 Cover Temperature, Run ID, Notes などをタッチすると、その他の情報 を入力することができます。 デフォルトの Run ID は年月日時間(YYYYMMDDHHMM)の順番で 表示されます。YYYYMMDD は日付で HHMM が現在時刻です。

- 6. ランを開始します。
  - 1台の単独の装置としてランを実行する Start Run Now をタッ チして、確認画面が現れたら OK を選択します。装置のヒーターカ バーが設定温度に達したらランが開始されます。ランが終了すると、 ランレポートが自動的に表示されます。
  - 複数の装置を同時にラン実行する Start Multiple Runs をタッチ すると、ネットワーク接続をしている装置が、Select Instrument 画面が表示されます。

Select Instruments	×
[Ready] [Ab2 [Ready]	lab3 [Incompatible]
Instrument1 (D_001 [Ready] [Disallowed]	
(Running) ID_002	
Instrument2 ID_003 [Ready]	
Number of hubs selected: 2	Start Run Now
Select the instruments for the run. To run the method on the current instrument, touch the button in the upper left corner.	?

- 7. Select Instrument 画面では
  - a. ランを実行する装置のチェックボックスにタッチしてチェック を入れます。この画面には実行するランメソッドが利用可能な 装置のみ選択可能で、利用ができない装置はチェックボックス が ⊗ で表示されます。 ランを実行したい場合は、画面の左上のチェックボックス (LocalHab)を選択します。
  - b. Start Run Now をタッチし、確認画面が現れますので OK を選択 します。ランメソッドが選択した各々の装置に送られ、装置に 問題が無い場合はランが実行されます。 ラン開始後、ランメソッドの変更、ポーズおよびランの停止を 行いたい場合は、リモートで運転している装置を直接操作する 必要があります。ランが終了すると、ランレポートが表示され ます。

## 2.6 ランのモニタリング

Run Monitor 画面 以下に示す Run Monitor 画面でランの状態を確認することができます



Run Monitor 画面の情報

	図示番号	記述
А	Run Monitor ボタン	現在のサンプル温度、ランの終了予定時間、
	とステイタスバー	現在実行中のステージを表示しています。
В	ホットアイコン	ヒーターカバーが火傷の恐れがある温度に
		達していることを示しています。
С	ステージ情報	VeriFlex <sup>TM</sup> システムが現在実行しているス
		テージとサイクル番号を示しています。
D	ステップ温度	ステップの温度を示しています。
Е	保持時間	ステップの保持時間を示しています。実行中
		のステップの場合は時間がカウントダウン
		されます。
F	ステップ情報	VeriFlex <sup>™</sup> システムが実行しているステッ
		プを示しています。
G	ステータスレポート	ランのステイタスやラン中に起きたエラー
		等を表示しています。
		右端の三角形をタッチすると、エラーレポー
		トの全文が表示できます。
Н	矢印ボタン	ランメソッドの全ステップが画面サイズで
		表示できなかった時に表示されます。矢印ボ
		タンにタッチするとランメソッドがスライ
		ドして、未表示部分を表示させることができ
		ます。

p41 の"ラン中の Run Monitor 画面における操作"を参照して、ラン中にで きる操作を確認してください。

### ラン中の Run Monitor 画面 における操作

ラン実行中でもタッチ画面を利用して表1の操作をすることができます

ラン中の Run Monitor	・画面における操作
------------------	-----------

操作	使用するキー
表示できていないステッ プを表示する	左右の矢印ボタンをタッチして画面をスク ロールします
ランの一時停止	<b>Pause Run</b> をタッチします("ランのポーズ" p3-19 を参照してください)
ランメソッド情報の閲覧	View Method Information をタッチします。入力したランのパラメーター情報が表示されます
ラン時間の表示	<ul><li>Display Method Time をタッチします。ランの経過した時間と残りの時間を表示します</li></ul>
次のステップにスキップ する	<ul> <li>Skip をタッチします。装置が温度の保持 を行っているときのみスキップ機能が利 用できます</li> </ul>
ランを停止する	Stop をタッチします。
サイクルステージのサイ クル数の変更	サイクルフィールドをタッチして、サイクル数の 数値を変更します("サイクル数の変更" p3-18を 参照してください)
Main Menu 画面を表示 する	ホームボタンをタッチします。ランの実行 中でも、Main Menu 画面から Veriti™ サ ーマルサイクラーのほとんどの機能にア クセスすることができます
ステータスレポートを表 示する	▲ ステイタスボタンをタッチします

- **サイクル数の変更** ラン実行中であってもサイクルステージのサイクル数を変更することができ ます。
  - 1. オレンジの Run Monitor ボタンで Run Monitor 画面に戻ります。
  - 2. サイクルステージのサイクルフィールドをタッチして、新しい値を入力 します。ただし以下の条件では変更はできません。
    - a. すでに完了しているステージの変更
    - b. すでに完了しているサイクル数よりも小さい値の入力
  - 3. Send をタッチすると Run Monitor 画面に戻ります。 新しく設定されたサイクル数でランは継続されます。

1. Run Monitor 画面で Pause Run をタッチします。



**p53**の "ポーズまたはスタンバイの持続時間"で設定された時間を最大 として、ランを一時停止することができます。 Pause Run ダイアログボックスでは2つの情報が表示されます。

- 一時停止の残り時間が Remaining Time として表示されます。
- 一時停止をしてから経過した時間が Elapsed Time として表示され ます。
- 2. 下記の操作をするには各々の操作を選択してください。
  - デフォルトで設定したポーズ時間を満了させる場合: 何もする必要はありません。
  - 特定の時間を設定して一時停止したい場合: Edit Time をタッチし、 一時停止の時間を入力します。そして Done を選択してください。 一時停止時間を満了するとランに復帰します。
  - すぐにランを復帰させたい場合: ダイアログボックスの左にある Resume Run をタッチします。
- **ランの停止** Run Monitor 画面で Stop とタッチした後、確認画面が現れますので OK を選 択してください。ランが終了して、ランレポートが表示されます。

Run Monitor ボタンを使う Run Monitor ボタンは Run Monitor 画面の上部に表示されています。もしラン中に Run Monitor 以外の画面へ表示を切り換えた場合でも、このボタンはすべてのメニュー画面の上部に表示されます。 このボタンをタッチすれば、どの画面からでも Run Monitor 画面に戻れます。

Sample:<br/>61.0 °CStage 1Time Remaining:<br/>00:01:17Run Monitor ボタン

もしラン中にエラーが発生した場合、ボタンの中央にある実行中のステージの 表示がアラートアイコンに変わります。



## 2.7 ランレポートの閲覧と印刷

ラン終了後、以下に示すランレポート画面が表示されます。もしランメソッド の最後のステージが無限大保存の設定の場合は、まず Stop ボタンをタッチし た後、確認画面で OK を選択するとランレポートが表示されます。 ツールメニューから View Last Run をタッチしても、ランレポートを表示する ことができます。

Run Re	port	×
	Date & Time: 2007-02-14 10:07:11 PM Instrument Name: 123456769012345676901234 Initiating Hub: 123456789012345678901234 Run Method: LMS2 Run ID: 200702140936 User Name: gregor Reaction Volume(uL): 10 Cover Temperature(Degree C): 105.0	
	Run Started: 2007-02-14 09:37:35 PM Run Completed: 2007-02-14 10:07:11 PM Notes: This is a quick test for Joe's samples with a 5 minute vortex step. Run Status:	Save

NOTE: ランレポートは次のランが完了するまで保存されます。

- ランレポートを印刷するには、ランレポート画面から Print をタッチします。もしいくつかのエラーがラン中に発生した場合は、エラー情報も同時に印刷されます。
- ランレポートをUSB ドライブに保存するには、Save をタッチします。

## 2.8 装置からサンプルを取り出す

サンプルトレー/リテーナーアセンブリ、または 96 ウェルプレートを両手でつ かみ、ゆっくり前後にゆすってブロックから外します。

!注意! サンプルブロックの温度が 27℃以上のときは、サンプルキャッ プがはじけるように開いてしまうことがありますのでご注意ください。

## 2.9 サイクルパフォーマンステストの実行

<< 目的 >>

このテストではサンプルブロックが適切に加熱・冷却されるかを検定するのと 同時に、PCR サイクルが適切に機能するかを検定します。

**必要なもの** テストを行なう際には、サンプルブロックに適合するプレートが必要です。

- Veriti<sup>™</sup> Fast サーマルサイクラーの場合: MicroAmp<sup>™</sup> Fast 96-Well Reaction Plate(商品番号 4346907)
- Veriti<sup>™</sup> サーマルサイクラーの場合: MicroAmp<sup>™</sup> 96-Well Reaction Plate (商品番号 N8010560)
- 操作手順 1. Main Menu 画面から Tools Menu を選択します。
  - 2. Tools Menu 画面で、Run Cycle performance Test を選択します。
  - 3. もしこのテストの結果を保存したい場合は、この時点で USB メモリを 挿入してください。
  - 4. 画面の指示にしたがって、テストを開始してください。テスト結果は "Pass" または "Fail" で表示されます。
  - 5. テストの結果が Fail の場合は、ヒーターカバーが閉まっていることを 確認して、もう一度テストをやり直してください。
  - 6. もし再度 Fail となった場合は、弊社および弊社代理店までご連絡くだ さい。
  - 7. テスト結果が Pass となれば、サイクルパフォーマンステスト OK です。

NOTE: このサイクルパフォーマンステストは約10分で終了します。

## 3. Veriti<sup>™</sup>サーマルサイクラーの初期設定

次の項目のデフォルト設定、またはセキュリティ機能を設定することができます。

- 日付と時間の設定 (p46)
- セキュリティおよびユーザーアカウントの設定 (p47)
- ヒートカバー及びサンプルブロックのアイドリング温度 (p52)
- ポーズまたはスタンバイの持続時間 (p53)
- Shortcut の設定: Main Menu 上に表示 (p53)

**日付と時間の設定** 日付と時間はランレポートに記載されるほか、装置の多くの画面の下部に表示 されます。

- 1. Main Menu 画面の Settings Menu をタッチします。
- 2. Settings Menu 画面の Admin Menu をタッチします。
- 3. Admin Menu 画面の Set Date&Time をタッチします。
- 日付をセットする場合は Date のフィールドを選択し、YYYY-MM-DD (年・月・日)のフォーマットで今日の日付を入力します。"- (ハイ フン)"は自動的に入力されます。
- 時間をセットする場合は Time のフィールドを選択し、HH:MM:SS (時:分:秒)のフォーマットで現在時刻を入力し、AM または PM を 選択します。":(コロン)"は自動的に入力されます。
- 6. **Done** を押した後、**OK** をすると設定条件が保存され、Admin Menu 画 面に戻ります。

## セキュリティおよびユーザーアカウントの設定

**セキュリティ設定** Veriti<sup>™</sup> サーマルサイクラーのすべてのセキュリティ機能は、デフォルトでは の有効化 オフになっています。この状態では、すべての人が装置を使用すること、およ び装置の各設定を変更することが可能です。

《装置のセキュリティ設定を変更》

- 1. ログインが必要な設定になっている場合は、Administrator 権限のユー ザーでログインします。
- 2. Main Menu 画面の Settings Menu をタッチします。
- 3. Settings Menu 画面の Admin Menu をタッチします。
- 4. Admin Menu 画面の Security をタッチします。

Security			×
Require Login:	Ves	No No	
Enable Guest Account:	Ves	No No	
Allow Empty Password:	Yes	V No	
		Done	
Touch the buttons to ena and/or a guest account, to accounts with no passwol	ble or disable login o allow/disallow user rd, then touch Done.		?

- 5. Security 画面で、各ユーザーが装置使用時にログインが必要にするに は、"Require Login"の Yes を選択します。
- ゲストがログインすることを許可する場合は、"Enable Guest Account"のYes を選択します。
   Guest ユーザーは、ランメソッドの実行・閲覧・プリントと各テスト の実行が可能です。但し、ランメソッドがUSBドライブで保存されて いる場合は、ランメソッドの作成・閲覧・編集・削除・保存・プリント が可能になります。
   ゲストユーザーは、ユーザーアカウントや各プリファレンス設定を管理 することはできません。
- 7. パスワードなし(ブランク)での新規ユーザーアカウントを許可する 場合は、"Allow Empty Password"の Yes を選択します。

Note: このオプションを Yes とした場合も、既にパスワードが設定されているユーザーについて は、ログイン時にパスワード入力が必要です。

- 8. Done をタッチした後、
  - Yes を選択すると、装置が再起動し、設定した変更点がすぐに有効になります(但し、ランが実行中の場合は、再起動は行なえません)。
     再起動後の操作については、p48の"新規ユーザーの設定"をご参照ください。
  - Noを選択すると、装置背面の電源スイッチによって次回再起動した時に変更内容が有効になります。
- **アクセスレベル** 各ユーザーは、装置をどこまでコントロールできるかを決定するアクセスレベ ルが割当てられています。アクセスレベルは次の3つに分かれます。
  - Limited User: ランメソッドの実行・閲覧・プリントと各テストの 実行が可能です。ランメソッドの作成・編集は行なえません。また、 ユーザーアカウントや各プリファレンス設定を管理することはできま せん。
  - User: ランメソッドの実行・作成・閲覧・編集・削除・保存・プリントと各テストの実行が可能です。ユーザーアカウントや各プリファレンス設定を管理することはできません。
  - Administrator: セキュリティ設定、ユーザーアカウントや各プリフ アレンス設定の管理を含め、装置の持つすべての機能にアクセスでき ます。
- 新規ユーザーの 装置がユーザーによるログインが必要な設定になっている場合、デフォルトの
   設定 Administrator ユーザーの名前とパスワードを使用してログインし、新たな利用者固有のユーザーアカウントを作成する必要があります。

《新規ユーザーの設定方法》

1. Veriti<sup>™</sup>サーマルサイクラーの起動画面で、On をタッチし、引続き Log In をタッチします。



- 2. 次の手順でログインします。
  - a. User フィールドに、"admin" と入力します。
  - b. Password フィールドに、"password" と入力します。
  - c. **Log In** をタッチします。
- 3. Main Menu 画面の Settings Menu をタッチします。
- 4. Settings Menu 画面の Admin Menu をタッチします。
- 5. ユーザーの設定については、次の"ユーザーの追加"を参照して下さい。 管理者としてはじめてユーザー設定を行なう場合は、必ず Administrator として作成して下さい。
- **ユーザーの追加** 《ユーザー設定方法》
  - 1. Administrator としてのアクセス権限でログインします。
  - 2. Main Menu 画面の Settings Menu をタッチします。
  - 3. Settings Menu 画面の Admin Menu をタッチします。
  - 4. Admin Menu 画面の Manage Users をタッチすると、Manage Users 画 面が表示されます。

Manage Users			×
User	Last Login	Access	
admin	2006-12-07	Administrator	
myuser	2006-12-07	User	
myadmin	2006-12-07	Administrator	Page 1 / 1
Add User Edit User Delete User			
Touch a user name to select it, then touch any of the buttons to perform an action. Touch a column title to sort the table.			?

5. Add User をタッチします。

Create New User Account 画面が表示されます。

User Name:*		
assword*:		
Re-enter Password:	*	
Phone:		
Email:		
Permission Level:*		
ast Update:		
	Set Access	Done

- 6. 新しいユーザーに関する情報を入力します。入力必須フィールドは、 User Name, Password, Re-enter Password です。それ以外のフィール ドはオプションです。
  - a. User Name は 5~15 文字で入力します。スペースは使用でき ません。
  - b. Password は、少なくとも1文字以上で設定します。Security 画面で "Allow Empty Password"にチェックが入っている場 合を除いては、スペースを入力するはできません。"Allow Empty Password"が有効な場合は、パスワードなし(ブラン ク)での設定が可能です。
  - c. Set Access をタッチし、次のユーザータイプから1つを選択し ます。
    - Limited User
    - User
    - Administrator
  - d. Done をタッチして Permission 画面を閉じます。
- 7. Done に続き、OK を押すと装置にユーザーアカウントが追加されます。 新しいユーザーが Manage Users 画面に現れます。
- 8. ユーザー設定が完了したら、× を押して、Admin Menu 画面に戻り ます。

ユーザーの 《ユーザーの変更・編集》

変更·編集

- 1. Administrator のアクセス権限でログインします。
- 2. Main Menu 画面の Settings Menu をタッチします。
- 3. Settings Menu 画面の Admin Menu をタッチします。
- 4. Admin Menu 画面の Manage Users を押すと、Manage Users 画面が表示されます。



5. 編集したいユーザーをリストから選択し、Edit User をタッチします。 Edit User Account 画面が開きます。

Edit User Acc	count 🛛 🗙
User Name:*	admin
Password*:	*****
Re-enter Password:*	****
Phone:	123456
Email:	tcadmin@appliedbiosystems.com
Permission Level:*	Administrator
Last Update:	
	Set Access Set Done Done
Update by touc	Account information accordingly ching on the relevant information

- 6. Edit User Account 画面上で、変更内容を入力します。
  - Phone または Email の入力フィールドを選択し、情報を入力します。
  - Set Access をタッチして、ユーザーのアクセスレベルを変更します。
  - Set Password をタッチして、ユーザーのパスワードを変更します。
- 7. Done に続き、OK をタッチすると Manage Users 画面に戻ります。
- 8. 設定が完了したら、× をタッチして、Admin Menu 画面に戻ります。
- **ユーザーの削除** 《ユーザーの削除方法》
  - 1. Administratorのアクセス権限でログインします。
  - 2. Main Menu 画面の Settings Menu をタッチします。
  - 3. Settings Menu 画面の Admin Menu をタッチします。
  - 4. Admin Menu 画面の Manage Users を押すと、Manage Users 画面が表示されます。



- 5. 削除したいユーザーをリストから選択し、Delete User をタッチします。 確認画面が表示されます。
- 6. **OK**を押すと、ユーザー、およびユーザーが作成したすべてのランメソ ッドとフォルダが削除されます。引続き、**OK**で決定します。
- 7. 設定が完了したら、× を押して、Admin Menu 画面に戻ります。

## その他のオプション設定

下表は調整または作動可能なオプション設定です。これらのオプションはすべてのユーザーで設定可能です。

Veriti™サーマルサイクラーのオプション設定と各デフォルト値

オプション	デフォルト値	ページ
ヒートカバー及びサンプルブロック	ヒートカバー: 105℃	52
のアイドリング温度	サンプルブロック:25℃	
ポーズまたはスタンバイの持続時間	ポーズ時間: 10 秒	53
	スタンバイ: 1 分 30 秒	
Main Menu 画面の"Shortcut"の設	なし	53
定		

- ヒートカバー及び
- 1. Main Menu 画面の Settings Menu をタッチします。

サンプルブロック のアイドリング 温度の設定

2. Settings Menu 画面の Set Idle Temperature をタッチします。



3. Idle State Temperature 画面で、ヒートカバー、またはサンプルブロックのアイドリング条件を設定します。

- a. Enabled を選択するとオプション設定が有効になります。
   Disabled が選択されていると、ヒートカバー、およびサンプル ブロックのアイドリング温度は装置によって維持管理されず、
   室温下での温度となります。
- b. 各 Temperature フィールドに設定温度を入力します。サンプル ブロック温度は 4~99℃の間で設定します。
- 4. Done を押すと変更内容が保存され、 Settings Menu 画面に戻ります。
- ポーズまたは スタンバイの 1. Main Menu 画面の Settings Menu をタッチします。

スノンハーの 持続時間

- 2. Settings Menu 画面の Set Time-outs をタッチします。
  - 3. Set Time-Out 画面上で、Time-outs の入力フィールドを選択し、持続 時間を入力します。
    - Default Standby Time-Out は、装置のアイドリング状態においてタ ッチ画面画面が消えるまでの持続時間を設定します。タッチ画面画 面が消えないようにするには、00:00 と入力します。
    - Default Pause Duration は、ラン実行中にポーズをかけた時のポーズの持続時間を設定します。
  - 4. **Done** をタッチした後、**OK** をすると設定条件が保存され、Settings Menu 画面に戻ります。

 Main Menu 画面の Shortcut の設定
 Main Menu 画面にある Shortcut ボタンを利用して、装置の様々な機能をダイ レクトに実行できるよう設定することが可能です。装置使用時にログインが必 要な設定の下では、各ユーザー毎に独自のショートカットを設定することがで きます。ログインが不要な場合は、すべてのユーザーに対して共通の Shortcut ボタンが表示されます。

- 1. Main Menu 画面から、次のどちらかのボタンをタッチします。
  - Shortcut X (X=番号)
  - Settings Menu  $\rightarrow$  Configure Shortcuts

2. Configure Shortcuts 画面にショートカットとして適用可能な機能リストが表示されますので、設定する機能を選びます。 ▲ と ◆ をタッチして、リストをスクロールします。

Configure Shortcuts	×
About the Instrument	
Set Default Method	Page 1/3
Set Instrument Name Number of selected shortcuts: 2	Done
Select the functions that will be shown on the main r	nenu. <b>?</b>

3. Done をタッチして、Settings Menu 画面に戻ります。 設定したショートカットが Main Menu 画面に表示されます。

### www.appliedbiosystems.co.jp

### ■弊社製品に関わるアプリケーション等の情報、技術資料ライブラリ

下記弊社ウェブサイトをご覧ください。 http:www.appliedbiosystems.co.jp

よくあるご質問: トップページ→Support→「製品Q&A」 技術資料ライブラリ: トップページ→Support→「テクニカルアップデート」

TEL. 0120-477-392 FAX. 0120-477-120

# ■弊社機器修理・保守契約等のご質問とご依頼TEL. 0120-203-885 FAX. 03-5566-6503

Applera Corporationは、ライフサイエンスに関する世界有数の技術と情報を提供しています。 Applera Corporationは、Applied BiosystemsとCelera Genomicsから構成されています。 © 2007 Applied Biosystems Japan Ltd., All rights reserved. Printed in Japan. 商品番号: JP 777



アプライドバイオシステムズジャパン株式会社

取扱店:

本社: 〒104-0032 東京都中央区八丁堀4-5-4 TEL.03 (5566) 6100 FAX.03 (5566) 6501

大阪: 〒564-0052 大阪府吹田市広芝町10-28 TEL.06 (6389) 1201 FAX.06 (6389) 1206