applied biosystems

VeritiPro[™] Thermal Cycler INSTALLATION, USE, AND MAINTENANCE

For 96-well and 384-well VeritiPro[™] Thermal Cyclers

for use with Invitrogen[™] and Applied Biosystems[™] PCR reagents

Catalog Numbers A48141, A48140

Publication Number MAN0019157

Revision B.0







Life Technologies Holdings Pte Ltd | Block 33 | Marsiling Industrial Estate Road 3 | #07-06, Singapore 739256 For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. MAN0019157

Revision	Date	Description
B.0	7 April 2021	Addition of specifications for 384-well model instrument.
A.0	11 September 2020	New document.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

TRADEMARKS: All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

©2021 Thermo Fisher Scientific Inc. All rights reserved.

Contents

	CHAPTER 1 Product information	7
	Product description	7
	Instrument overview	7
	CHAPTER 2 Start, sign on, and configure the instrument	. 8
	Required materials not provided	8
	Set up the VeritiPro [™] Thermal Cycler	8
	Touchscreen controls	. 11
	Home screen controls	. 12
	Enter text	. 13
	Enter numbers	. 13
•	CHAPTER 3 Connect the instrument for remote monitoring (Optional) Set up the wired connection	. 14
	Create a user profile on the PCR instrument	. 18
	Create a Connect account	
	Link the PCR instrument to Connect	
	Connect by mobile device	
	Connect by PC	
	Connect by instrument	. 20
	Manage user profiles	. 21
	CHAPTER 4 Methods	22
	Create/Edit Method	. 22
	Create a new Method	
	Edit a method	
	Save a Method	. 25

Actions	26
Manage Methods and Folders	32
Open and run Method	. 36
Load samples into the instrument	36
Perform a run	. 36
Monitoring a run	. 39
View and export the run report	. 41
Remove the samples from the instrument	41
CHAPTER 5 Maintenance	. 42
Clean the instrument	. 42
Preparation	
Clean the touchscreen	
Clean the sample wells	. 43
Decontaminate the sample wells	43
Clean the heated cover	43
Decontaminate the heated cover	. 44
Replace the fuses	45
Required materials	45
Procedure	. 45
Upgrade the system firmware	46
Determine firmware version on instrument	. 46
Upgrade the instrument firmware (Cloud)	. 46
Download new firmware	. 47
Upgrade the instrument firmware (USB drive)	. 47
Self Verification test	. 48
Procedure	. 48
Restore factory settings (Administrator profile only)	49
APPENDIX A Troubleshooting	. 50
Troubleshooting	50
Return an instrument for service	. 52
APPENDIX B Recommended instrument settings	53
About Instrument	. 53
Recommended instrument settings	. 53
Instrument settings	
Maintenance and services	
Run history	
LISB shortcuts	60 60

	APPENDIX C	Instrument specification and layout	61
	Site rec	quirements	61
	Environ	nmental requirements	61
		emperature and humidity requirements	
		bllution	
		titude	
	•	cations	
		echnical specifications	
		and communication port symbols	
		on of power point and ports on the instrument	
	Locatio	on or power point and ports on the instrument	04
	APPENDIX D	Predefined Methods	65
	Predefi	ined run methods	65
-	APPENDIX E	About the Thermal Cycler Fleet Control Software	88
	Connec	ct the thermal cycler to the network	88
	APPENDIX F	Ordering information	89
	Instrum	nent part numbers	89
	Consur	mables	89
	APPENDIX G	Safety	92
	Symbo	ols on the instrument	92
	Safety	symbols on this instrument	93
	Lo	ocation of safety labels on the instrument	94
	China F	RoHS Compliance	95
	Safety	information for instruments not manufactured by Thermo Fisher Scientific	96
	Instrum	nent safety	96
		eneral	
		nysical injury	
		ectrical safety	
		eaning and decontamination	
	•	and electromagnetic compatibility (EMC) standards	
		afety	
		ass A notice	
		nvironmental design standards	

Chemical safety	99
Biological hazard safety 1	100
APPENDIX H Documentation and support	01
Related documentation1	101
Customer and technical support1	101
Limited product warranty 1	101



Product information

Product description

The VeritiPro[™] Thermal Cycler is an end-point thermal cycler, specifically designed for the amplification of nucleic acids using the Polymerase Chain Reaction (PCR) process. The user interface includes a touchscreen with a graphical display that shows the time, status, and temperature for each run. A touchscreen keypad allows you to enter information into fields on the display screen. It is also compatible with the Applied Biosystems[™] Thermal Cycler Fleet Control Software.

There are two different models of the thermal cycler, one with a 96-well VeriFlex[™] block (Cat. No. A48141), and the other with a 384-well block (Cat. No. A48140).

Instrument overview

The VeritiPro[™] Thermal Cycler allows you to:

- Optimize your PCR with the help of the VeriFlex[™] block (96-well block only)
- Program the instrument from the touchscreen interface
- · Access the system remotely through a mobile application
- Simulate other PCR instruments with Thermal Simulation modes



Figure 1 VeritiPro[™] Thermal Cycler

This user guide provides information on how to maximize the capabilities of this system.

For detailed specifications of the VeritiPro[™] Thermal Cycler, see Appendix C, "Instrument specification and layout".



Start, sign on, and configure the instrument

Required materials not provided

- Can of compressed air (For use in cleaning wells of sample block).
- (Optional) Electrical protective devices.

Note: The use of one or more of the following electrical protective devices is recommended.

- Power line regulator (100-240 V)
- Surge protector/line conditioner (10-kVA)
- Uninterruptible power supply (1.5-kVA)

Set up the VeritiPro[™] **Thermal Cycler**

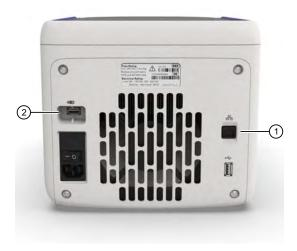
IMPORTANT! Save the packing materials and box in case you need to ship the instrument in for service.

- **1.** Open the shipping crate to unpack the instrument. You should receive one box containing the thermal cycler and the accessories.
- 2. Remove the packing material, then inspect the instrument for shipping damage.
- **3.** Use compressed air to clear out each well of the sample block to remove particles that may have collected inside during shipping.
- **4.** Move the instrument to an installation site that meets the spatial and weight requirements for the thermal cycler (see "Specifications" on page 62).

- **5.** Connect the instrument to a power outlet.
 - a. Connect the power cord to the instrument.



- 1 Power cord port
- **b.** Install any desired electrical protective devices.
- c. Connect the power cord to a wall plug.
- **6.** Connect the instrument to a network.
 - a. Connect the ethernet cable to the ethernet port on the instrument.
 - b. Connect the ethernet cable to the ethernet port wall plug.
 In the absence of the ethernet cable or the ethernet port, connect a High-Power USB Wi-Fi Module (Cat. No. A26774) to the instrument using the USB port on the instrument.



- 1 Ethernet port
- 2 USB port for wi-fi module



7. Press the power switch, then wait for the instrument to start up. Proceed with the installation after the touchscreen displays the Home screen, indicating that the instrument is active.



1 Power switch

When you power on the instrument, the instrument will take under a minute to start up. The Home screen will display following successful installation.



Touchscreen controls

Table 1 General touchscreen controls

Button	Function
•	Returns to the previous screen
	Go to Home screen
	Go to Sign in screen
*	Go to Settings screen
> <	Advance and return through stages
×	Close the current modal window.



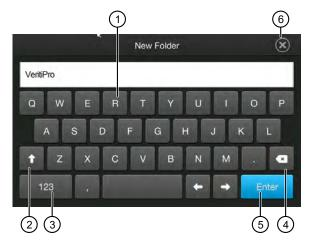
Home screen controls

Table 2 Home screen controls and indicators

Button	Function
	Go to Set up run screen
45.2°C ①	View instrument status
01:02:23	1. Current block temperature
Running 3	2. Time remaining
A CONTRACTOR OF THE PARTY OF TH	3. Status of the block (pre-heating, running, etc.)
	Go to Sign in screen
	Go to Settings screen
<u></u>	Indicates whether the Wi-Fi is on or off.
	Indicates whether a USB device is inserted into the instrument.
윰	Indicates whether the instrument is connected to the network.

Enter text

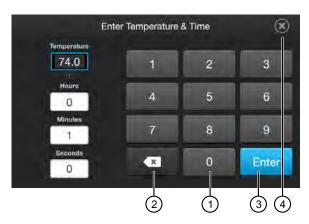
When you press a field that requires the input of text, the text editor, as seen in the following figure, opens.



- 1 Enter a letter
- (2) Change letter case
- (3) Enter punctuation or other symbols
- 4 Delete
- (5) Close and save
- 6 Close without saving

Enter numbers

When you press a field that requires a numerical input, the numeric editor, as seen in the following figure, opens.



- 1 Enter a number
- 2 Delete or enter decimal, colon or infinity
- (3) Close and save
- 4 Close without saving



Connect the instrument for remote monitoring (Optional)

Set up the wired connection

See Chapter 2, "Start, sign on, and configure the instrument" for details on connecting the instrument to the network. To set up the wired connection:

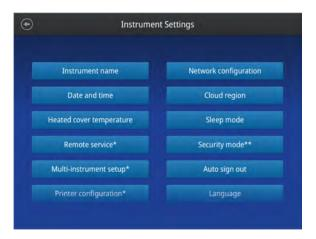
1. On the **Home** screen, select **(Settings)**.



2. In the Settings screen, select Instrument Settings.



3. In the Instrument Settings screen, select Network configuration.



4. In the **Network Connection** screen, select a field in the **Wired** panel.



- (1) Wireless panel
- (2) Wired panel
- 5. Select a method to enter an IP address.
 - **a.** Select **DHCP** to obtain an IP address automatically. A check mark appears when DHCP is selected.



b. Select **Static IP** to enter an IP address manually, then enter the appropriate IP addresses for the instrument, the Subnet Mask, and, optionally, the Default Gateway, the Primary DNS Server, and the Secondary DNS Server using the numeric editor. Addresses are in the form of X.X.X.X, where each X is a 3- digit number, from 001 to 255.



Note: If your instrument is not on a network, you do not need to set the IP address. Ask your system administrator if the IP address is assigned statically or dynamically. For static addresses, you need to know the IP address for the instrument, the subnet mask, and the default gateway.

6. Select **Done** to save the changes and go back to the **Instrument Settings** screen or select **Cancel** to exit the screen without saving the changes.

Set up the wireless connection

See Chapter 2, "Start, sign on, and configure the instrument" for details on connecting the High-Power USB Wi-Fi Module to the instrument.

- **1.** See "Set up the wired connection" on page 14 Steps 1 through 3 to find the **Network configuration** screen.
- 2. In the Network configuration screen, select a field in the Wireless panel.



- (1) Wireless panel
- (2) Wired panel

Note: During initial setup, if you selected the Wired option in the **Network Connection** screen, you will be required to enter the IP address if you selected the Static IP wired option. If you selected the Dynamic IP wired option, the IP address is automatically populated.

3. Once a wireless connection has been detected, a list of the available networks is displayed. Select the network name of your choice or select **Join others**.

Note: If you choose Join others, the Configure and Join Network screen opens.

- **4.** In the **Configure and Join Network** screen, select the **Network Name** field, then enter the name and security type of the network.
- **5.** Select the security type from the **Security type** drop-down menu.

Note: Contact your IT Systems Administrator for information on security type.

Select from the following options:

- Open
- WEP
- WPA Personal
- WPA2 Personal

- WPA Enterprise
- WPA2 Enterprise

Note: The above options are available only if Join Other Network was selected in Step 3. You cannot change the security type if you selected an existing network.

- 6. Select Join to continue or Cancel to exit from the Find and Join a Network screen.
- 7. Depending on the security type you have selected, enter the appropriate passwords and select
- 8. If all the entered information is correct, the Network Connection Complete screen will appear. Select **OK** to continue.

Note: If incorrect information was entered the Network Connection Failed screen will open. Select **OK** to continue to the **Security type** screen.

Create a user profile on the PCR instrument

- **1.** Select . (Sign In) > Get started > Create profile.
- 2. Fill in the required text fields and enter a four digit PIN to create your user profile.

Note: The first profile created is automatically given an Administrator profile (indicated by an asterisk after the Username).

Create a Connect account

- **1.** Go to thermofisher.com/connect from your web browser.
- 2. Click **Sign up now** and follow the prompts to create an account.

Link the PCR instrument to Connect

- 1. Select . (Sign In) > Connect, then select the cloud region of the instrument.
- 2. Select the method for linking the instrument to Connect.

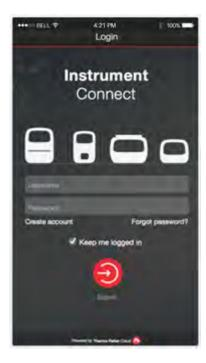


Connect by mobile device

Select **(Sign In)** • Get started • Connect • Mobile device from the instrument to generate a QR code.

- 1. Download the "Instrument Connect Mobile Application" on your mobile device.
 - **a.** For iPad[™] or iPhone[™] devices, download the application from the iTunes[™] music store by searching for Instrument Connect by Thermo Fisher Scientific.
 - **b.** For Android devices, download the application from Google[™] Play by searching for Instrument Connect by Thermo Fisher Scientific.





3. Capture the QR code on the instrument screen.

Connect by PC

Select & (Sign In) > Get started > Connect > PC from the instrument to generate a linking code.

- **1.** Log in to your Connect account using a web browser from a computer.
- **2.** Select (InstrumentConnect) from the left navigation strip.
- **3.** Select **(Add an Instrument)** from the top navigation strip.
- **4.** Select **VeritiPro** from the drop down menu, then click **Next**.
- **5.** Enter the linking code generated by the instrument in the text box, then click **Send**. Upon successful authentication, the instrument is linked to Connect.

Connect by instrument

- **1.** Enter your Connect **Username** and **Password** from the instrument.
- 2. Click Link account.

Note: If you do not have a PIN, you will be prompted to create one.

Upon successful authentication, the instrument is linked to Connect.

Manage user profiles

All users can manage their profiles to edit personal folder names, change PINs, and link to the cloud by selecting their **&** (Profile) to enter their My Profile page.

Users with Administrator profiles (as indicated by an asterisk after their user name) also have the ability to manage all user accounts by selecting **All accounts** after entering their **My Profile** page. Administrator have the ability to perform the following functions.

- · Create new user profiles.
- Grant Administrator privileges to user profiles.
- Manage all user folders.
- Delete PINs (this function effectively resets a PIN, so the user with a deleted PIN will need to create a new PIN the next time they log in).
- Delete user accounts.

Methods



Create/Edit Method

- Select any field on the graphical method profile to edit the parameter (temperature, time, heated cover, number of cycles, and volume).
- Select **Actions** (see "Actions" on page 26) to manage the steps, stages, and advanced options.
- Select Cancel to exit without making changes.
- Select **Save** as/**Save** to save new/updated method.
- Select Start Run to run method.

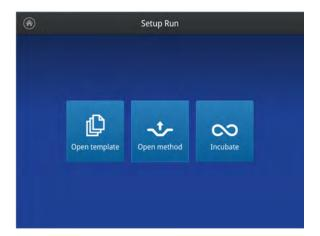


Create a new Method

1. In the Home screen, select Set Up Run.



2. In the Setup Run screen, select ((Open Template).



3. Select a folder , then select an appropriate template from the **Method Name** list.



4. Select **Actions** (see "Actions" on page 26) or any field on the graphical temperature profile to start editing the parameters.



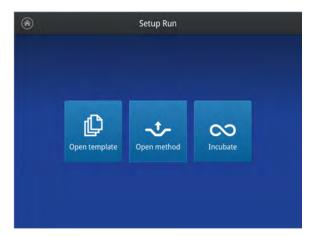
- 5. Select Save as, see "Save a Method" on page 25.
- 6. (Optional) Select Cancel to exit without making changes, or Start Run to run method.

Edit a method

1. In the Home screen, select Set Up Run.

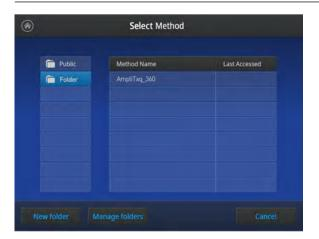


2. In the Setup Run screen, select - (Open Method).



3. Select a folder **a**, then select an appropriate template from the **Method Name** list.

Note: At least one method must have previously been set up in order for the list to be populated.



4. Select **Actions** (see "Actions" on page 26) or any field on the graphical temperature profile to start editing the parameters.



- 5. Select Save, see "Save a Method" on page 25.
- 6. (Optional) Select Cancel to exit without making changes, or Start Run to run method.

Save a Method

1. Once edits to a run method are complete, select **Save As/Save** to save the method.



2. In the Save As/Save screen, enter a name for the edited run method.

Characters allowed	Characters not allowed
<100 characters	>100 characters
Letters, numbers, spaces, underscores, and dashes	% * ? ; : , ! @ # \$. () < > / \ " ' ` ~ { } [] = + & ^ (tab)

3. Select **a**, then select the destination folder where you want to save the run method, then select **Save**.

Actions

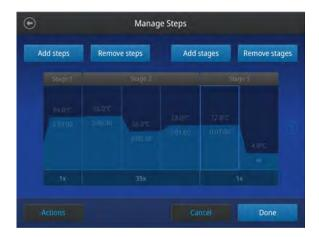
Select **Actions** to edit the parameters of a method. Options can be used to:

- Add or remove stages and steps
- Perform temperature optimization (manage VeriFlex[™] block zones, 96-well block only)
- Change ramp rates
- Increase or decrease time and temperature (manage AutoDelta)
- Run in simulation mode
- Add comments



Manage stages and steps

1. Select Stages and steps to open the Manage steps screen.



2. Select an option to manage stages and steps.

Press	Description
Add/Remove Steps, then - or +	To add/delete a step
Add/Remove Stages, then - or +	To add/delete a stage

- 3. Select **Done** when finished, or **Cancel** to exit the screen without making changes.
- **4.** Select **Save** (see "Save a Method" on page 25) to finish, or **Actions** to make additional changes.

Temperature optimization (96-well block only)

- **1.** Select **Temperature optimization** to edit the VeriFlex[™] block parameters.
- 2. Select r to choose the step to modify.



3. Select **Min/Max** to set a temperature gradient across the block, then enter values in the highlighted text fields.

Note: The maximum difference between the minimum and maximum temperature across the block is 30°C.



Chapter 4 Methods Create/Edit Method

4. Select **Midpoint** to set a midpoint temperature and temperature difference between zones, then enter values in the highlighted text fields.

Note: The maximum temperature increment or decrement cannot exceed 5°C.



5. Select **6-zones** to set the temperature for each zone of the block, then enter values in the highlighted text fields.



- 6. Select **Done** when finished, or **Cancel** to exit the screen without making changes.
- 7. Select Save (see "Save a Method" on page 25) to finish, or Actions to make additional changes.

Ramp rates

1. Select **Ramp Rates** to edit the ramp rate of the individual steps.



- 2. Select a ramp rate to open that ramp rate editor, then edit the ramp rate.
- **3.** Select **Done** when finished, or **Cancel** to exit the screen without making changes.

Note: If a simulation mode is set, then the ramp rate will be locked and unavailable for editing.

4. Select Save (see "Save a Method" on page 25) to finish, or Actions to make additional changes.

AutoDelta

- 1. Select Inc/Dec time and temperature to apply AutoDelta settings to a step.
- 2. Select to choose a step.



- 3. Enter the information for setting up AutoDelta.
 - **a.** Use the '-' and '+' toggle to set the Delta values temperature and time to be either decreased or increased.

Chapter 4 Methods Create/Edit Method

- **b.** Enter the cycle for the start of AutoDelta in the **Start AutoDelta from cycle** field.
- **c.** Enter the values for Delta temperature and/or time in the appropriate text fields.



- **4.** Select **Done** when finished, or **Cancel** to exit the screen without making changes.
- **5.** Select **Save** (see "Save a Method" on page 25) to finish, or **Actions** to make additional changes.

Remove AutoDelta

If AutoDelta has been applied to a step, a button will appear to allow removal.

Select Remove to disable AutoDelta.

Note: Temperature and time values will be cleared when AutoDelta is disabled.



Simulation mode

1. Select Simulation Mode to set the simulation mode.

Table 3 Simulation modes available on the VeritiPro™ Thermal Cycler

VeritiPro [™] 96-well Thermal Cycler	VeritiPro [™] 384-well Thermal Cycler
 No Simulation (default) Applied Biosystems[™] Veriti[™] Thermal Cycler Bio-Rad[™] T100[™] (Algo mode) Bio-Rad[™] MyCycler[™] BIOER[™] XP Cycler (Tube mode) Eppendorf[™] Mastercycler[™] MJ Research[™] PTC-200 Takara[™] Dice PCR Thermal Cycler Applied Biosystems[™] 2720 Thermal Cycler (Sample mode) GeneAmp[™] PCR System 9700 Bio-Rad[™] C1000 Touch[™] Biometra TAdvanced Eppendorf[™] X50a Intermediate GeneAmp[™] PCR System 9600 	 No Simulation (default) Applied Biosystems[™] Veriti[™] Thermal Cycler GeneAmp[™] PCR System 9700

Note: Simulation Mode makes this instrument perform like another instrument. This is done by simulating the ramp rate of another instrument. Not all of the listed simulation modes are available on all configurations of this instrument.

2. In the Simulation Mode screen, select the instrument you would like to simulate.



Note: Select **No Simulation** to run the instrument in the default mode.

3. Select **Done** when finished, or **Cancel** to exit the screen without making changes.

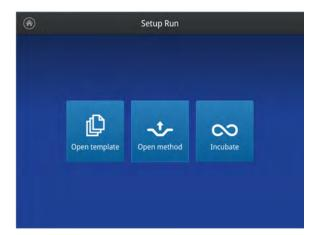
Manage Methods and Folders

Create a New Folder

1. In the instrument Home screen, select Set Up Run.



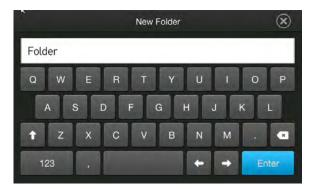
2. In the Setup Run screen, select - (Open Method).



3. In the Open Method screen, select New Folder.



4. Type a name for the new folder, then select **Enter**.



Manage Folders

If the instrument is linked to Connect, it will automatically display a Cloud folder.

IMPORTANT! Back up your method files to a USB drive frequently, or before sending in your instrument for service.

1. In the instrument Home screen, select Set Up Run → ❖ (Open Method) or ❖ (Open Method) directly from the bottom of the screen.



Chapter 4 Methods Create/Edit Method

2. In the Open Method screen, select Manage Folders.



3. In the **Manage Folders** screen, select folders from the left column and methods from the right column, then select one of the following options.

Select	Description	
Сору	To copy a run method from one existing folder to another	
Rename	To rename a particular method/ folder name	
Delete	To delete a particular method/ folder	



Copy a method

1. Select the **Method** to be duplicated.



- **2.** Select **Copy**, then paste the method to an existing folder or a new folder.
 - Select the destination folder for the duplicated method, then select **Paste method**.
 - Select New Folder and create a new folder for the duplicated method, then select Paste method.



3. Select Done when finished.

Open and run Method

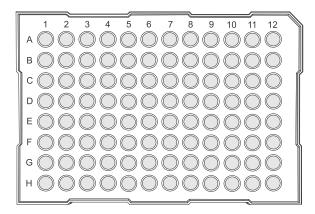
Load samples into the instrument

Note: Refer to the PCR Starter Kit user guides (see Appendix H, "Documentation and support") for instructions on filling, sealing, and loading MicroAmp[™] reaction plates, tube strips, and tubes.

Place the sample tray or plate onto the sample block(s)

The steps for placing the sample tray on the sample block(s) are the same for:

- A sample tray/retainer
- A sample tray without a retainer
- A 96-well or 384-well plate.
- **1.** (96-well block only) Lift the sample tray from the splash-free 96-well base and place it in the sample block(s).
- **2.** Place the MicroAmp[™] Tray or Plate onto the sample block(s) so that well A1 is at the upper left corner of the tray, as shown in the following figure:



3. Close the heated cover.

Perform a run

Prepare your samples and load the sample block(s) as described in "Load samples into the instrument" on page 36.

- 1. Close the cover.
- 2. Start the run from the **Home** screen in one of the following ways. Select:
 - (Open Template), see "Create a new Method" on page 22
 - (Open Method), see "Edit a method" on page 24
 - ° ∞ (Incubate)

3. Select Next, to display the Run Parameters screen.



4. (Optional) In the **Run Parameters** screen, select the field under **Verify the run ID**, then enter the applicable information in the **Enter the Run ID** screen.

Note: The default Run ID is YYYYMMDDHHMM, where "YYYYMMDD" is the date and "HHMM" is the current time.

- **5.** (Optional) Select **Comment** to enter any comments about the run.
- 6. Select Start Run.

The cover is heated to the required temperature, then the run begins.

While the run is in progress, the **Home** screen is displayed. When the run is complete, the **Status Dial** on the **Home** screen displays **Done** and **Remove Samples**.

7. Remove the samples and select **Done** after the run has finished.

Perform a run on multiple instruments

- 1. Select (*) (Settings) > Instrument Settings > Multi-Instrument Setup.
- 2. In the **Multi-Instrument Setup** screen, use the **Off/On** toggles to make the instrument visible and to find other instruments.

Option	Description
Make visible	Use the Off/On toggle to make the instrument on which you are running an experiment visible to other instruments on the network
Find other instruments	Use the Off/On toggle to find other instruments on the network.

- **3.** Select **△ ▶ △ ▶ ⑥** to return to the **Home** screen.
- **4.** Select **Set Up Run** and select **(Open Template)**) or **(Open Method)** and proceed to the **Start Run** screen. See "Perform a run" on page 36
- 5. In the Start Run screen, select Scan devices.

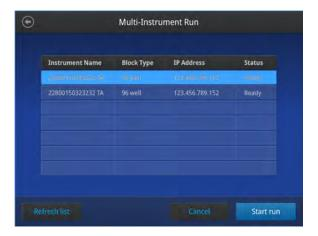
Note: The **Scan devices** button appears in the **Start Run** screen when you enable the multi-instrument run.



Note: The instrument starts scanning for all available VeritiPro[™] Thermal Cycler instruments on the same network which have the "Make Visible" setting enabled. All available instruments are displayed in a table in the **Multi-Instrument Run** screen.



6. Select one or more instruments from the table to enable the **Start Run** button.



7. Select Start Run.

Note: This machine cannot monitor runs taking place on other machines.

8. Remove the samples and select **Done** after the run has finished.

Monitoring a run

Monitoring a Run

• View the current run information on the **Home** screen dial according to the following.



- 1) Current block temperature
- (2) Time remaining since the run started
- 3 Run status
- Monitor the run progress in the Run Monitoring screen by pressing the Home screen dial.

Note: The **Run Monitoring** screen shows the run method profile and additional information that does not appear in the Dial on the **Home** screen.

Chapter 4 Methods Open and run Method



- 1) Start time displays the time a run started.
- 2 Time remaining displays the time remaining for a run.
- 3 End time displays the approximate time at which the run will end.
- 4 Details provides more details about the run.
 - Run ID
 - Run User
 - Simulation Mode
 - Start time
 - Estimated End time
 - Sample Temperature
 - Cover Temperature
 - Reaction Volume
- (5) Edit
 - Helps edit run parameters like temperature, time, and cycles.
 - Helps skip the current holding step. However, if the run is currently ramping, the Skip button will not function.
- 6 Pause to pause the run.
- 7 Stop run to stop the run.

Stop a run

- 1. Press the dial in the center of the screen.
- 2. Press Stop Run ➤ OK.
- 3. Remove the samples and press Done after the run has finished.

View and export the run report

- 1. Insert a USB memory device into the port on the front of the instrument.
- 2. In the Home screen, select **③** (Settings) ➤ Run History.
- 3. Select a run report to view.
- **4.** (Administrator profile only) Select **Manage ▶ Select all** to select all run reports. This option is not compatible with the print function.
- 5. Select Export to save the report on a USB memory device.
- **6.** (Optional) Select **Print** to print a single run report.

Note: You may be prompted to enter the IP address of the printer if it has not already been set.

Remove the samples from the instrument



CAUTION! During instrument operation, the temperature of the heated cover can be as high as 110°C, and the temperature of the sample block(s) can be as high as 100°C. Before performing the procedure, keep hands away until the heated cover and sample block(s) reach room temperature.



CAUTION! Sample caps may pop off if the cover is opened when the sample block(s) temperature is above 27°C.

Grasp the sample tray/retainer assembly or reaction plate firmly with both hands (or by using MicroAmp[™] Multi Removal Tool, Cat. No. 4313950) and gently rock it back and forth until the tubes (or plate) are free.

5

Maintenance

Clean the instrument



WARNING! Always wear protective glasses and gloves when servicing the instrument. Also, make sure you disconnect the instrument from AC line power before you begin any service procedure.

Preparation

Before cleaning the instrument:

- **1.** Power off the instrument by disconnecting the power.
- 2. Allow the instrument to cool until the heated cover and sample block(s) reach room temperature.

Clean the touchscreen

Clean the touchscreen with any commercially available LCD cleaning product. Be careful not to scratch the screen.

Clean the sample wells

If you use any cleaning or decontamination method, except those recommended in the manual, you risk damaging the equipment. Clean the sample wells once a month or as needed.



WARNING! During instrument operation, the temperature of the heated cover can be as high as 110°C, and the temperature of the sample block(s) can be as high as 100°C. Before performing the procedure, keep hands away until the heated cover and sample block(s) reach room temperature.

Clean the sample wells using a cotton swab dampened with isopropanol. **Do not** flush sample wells using a wash bottle or allow liquid to overflow from the wells.

- **1.** Follow the steps in "Preparation" on page 42.
- 2. Open the heated cover.
- 3. Remove the sample tray from the sample block(s) and set it aside.
- **4.** Use a cotton swab soaked in isopropanol to clean the sample wells thoroughly. Make certain that the isopropanol has evaporated completely before reloading a sample tray.

Decontaminate the sample wells

If the sample wells become contaminated with any biological agents, clean the wells thoroughly with a cotton swab soaked in 1:10 v/v dilution of 5.25% sodium hypochlorite (Clorox® bleach).

IMPORTANT! Use bleach solution in moderation. 10% bleach solution is recommended for removing contamination from the sample block(s), but excessive use of the solution can corrode the sample block(s) material. To prevent damage to the sample block(s):

- Avoid applying excessive amounts of bleach solution. If possible, instead of using a squeeze bottle or soaked cloth, use an atomizer to deliver the solution to the wells of the block(s).
- After treating with bleach solution, rinse the sample block(s) thoroughly using deionized water.

Note: Removing residual bleach from the surfaces of the instrument using water minimizes the long-term effects of bleach treatments.

If the sample block(s) become contaminated with radioactivity, use a commercially available decontaminant to remove the contamination. If the block(s) cannot be decontaminated, the instrument cannot be returned for service.

Clean the heated cover

Clean the heated cover once a month or as needed.

To clean the heated cover:

- **1.** Follow the steps in "Preparation" on page 42.
- 2. Open the heated cover.

Chapter 5 Maintenance Clean the instrument

3. Soak a cotton swab or piece of clean cloth with isopropanol and gently wipe the heated platen.



WARNING! CHEMICAL HAZARD. Isopropanol is a flammable liquid and vapor. Exposure may cause eye, skin, and upper respiratory tract irritation. Prolonged or repeated contact may dry skin and cause irritation. Exposure may cause central nervous system effects such as drowsiness, dizziness, and headache. Read the MSDS, and follow the handling instructions.

4. Remove any remaining isopropanol from the cover.

Note: If the platen becomes contaminated with amplified DNA, then raise the heated cover to the cleaning position, wipe the platen with a cloth or cotton swab soaked in bleach, then rinse with water.

Clean the heated platen once a month or as needed.

Decontaminate the heated cover

If the heated cover become contaminated with amplified DNA, raise the heated cover to the cleaning position and wipe the cover with a cloth or cotton swab soaked in 1:10 v/v dilution of 5.25% sodium hypochlorite (Clorox® bleach), then wipe the cover with a damp cloth.

Replace the fuses

Required materials

- Two UL Listed fuse, rated 10 A, Time-Lag T, 250 VAC, size: 5 × 20 mm
- Fine flat-tip screwdriver

Procedure



DANGER! ELECTRICAL SHOCK HAZARD. Severe electrical shock, which could cause physical injury or death, can result from working on an instrument when the high voltage power supply is operating. To avoid electrical shock, disconnect the power supply to the instrument, unplug the power cord, and wait at least 1 minute before working on the instrument.

- **1.** Power off the instrument by disconnecting the power.
- 2. Insert a small flat-tip screwdriver into the slot in the upper portion of the power entry module, and open the door.
- 3. Pull out the fuse compartment to inspect the two fuses in the fuse compartment:
- **4.** Pull out the blown fuse(s) from the fuse compartment and replace with new Two UL Listed fuse, rated 10A, 250VAC, size: 5 x 20 mm.
- **5.** Place the fuse compartment back into the power entry module (there is only one orientation) and close the door.
- **6.** Press the door until it locks in place.
- **7.** Connect the instrument power cord.

Upgrade the system firmware

Update software directly through Connect or using an USB drive with updated software downloaded from thermofisher.com/connect.



Determine firmware version on instrument

When a new firmware version is released, you may be required to load the new firmware on the instrument.

You will need a USB memory device and, if your instrument requires login, the login details to upgrade the firmware.

- 1. Select (3) (Settings) > About Instrument.
- 2. View current firmware version.

Upgrade the instrument firmware (Cloud)

IMPORTANT! You cannot upgrade the firmware while a run is in progress.

- 1. Select ② (Settings) > Maintenance & Services > Software Update > ThermoFisher Connect.
- 2. Select Yes to start the upgrade.

IMPORTANT! To prevent instrument malfunction and required service, do not power off the instrument during the upgrade.

When the upgrade process is complete, the instrument will automatically restart.

Download new firmware

- 1. Go to thermofisher.com from your web browser.
- 2. Click Product Support > Technical Resources > Product Support > Software, Patches & Updates.
- 3. Select VeritiPro[™] Thermal Cycler in the list, then click **Updates & Patches**.
- 4. Find the appropriate file. If the version number is:
 - The same as the current version on the instrument, you do not need to upgrade the firmware.
 - Different from the current version on the instrument, download the new firmware.
- **5.** Insert a USB memory device into the USB port on the computer.
- **6.** Click the link in the Software column, then select the USB memory device as the location for the saved file.

Note: The file must be downloaded to the root directory of the USB memory device and not into a folder.

7. Remove the USB memory device from the computer when the download is complete.

Upgrade the instrument firmware (USB drive)

IMPORTANT! You cannot upgrade the firmware while a run is in progress.

1. Insert the USB memory device with the new firmware in the USB port of your instrument.

Note: For instruments with the USB shortcuts feature enabled, you will be directed to the **USB** shortcuts screen. Select **Update Software** to proceed to the **Software Update** screen.

2. Select ② (Settings) ➤ Maintenance & Services ➤ Software Update ➤ USB drive.
The Software Update screen opens:



- 3. Choose the row with the new firmware file from the USB memory device, then Select.
- **4.** Select **Yes** to start the upgrade.

IMPORTANT! To prevent instrument malfunction and required service, do not power off the instrument during the upgrade.

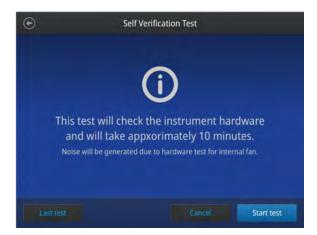
When the upgrade process is complete, the instrument will automatically restart.

Self Verification test

Use the **Self Verification Test** feature to check the instrument hardware. The check includes testing the block, heated cover, and other components.

Select Last Test to view the results of the last Self Verification Test.

Carry out the **Self Verification Test** periodically or whenever there is an intermittent instrument error. Contact your service representative in case of block failure.



Procedure

- 1. In the Home screen, select Settings > Maintenance & Services > Self Verification Test
- 2. In the Self Verification Test screen, select Start Test to begin testing or Cancel to exit the screen.

Note: The test takes about 10 minutes. Once the test is completed the test results will be displayed in the form of a report.

3. Select Export to save the test report to a USB drive or Close to exit the screen.

Restore factory settings (Administrator profile only)

Select **Restore factory settings** to remove all the data and customized settings and revert to factory settings. All data and settings will be erased once factory settings are restored. At the end of the restoration process, the message, "Your instrument has been restored." is displayed and the instrument automatically reboots after 30 seconds.





Troubleshooting

Troubleshooting

Use the following table to help troubleshoot problems with your Thermal Cycler.

Observation	Possible cause	Recommended action		
Instrument reports error(s).	Hardware malfunction.	Run the self verification test (see page 48).		
Reduction in instrument performance due to incorrect	Temperature sensor out of calibration.	Run a cycle performance test using the external multi-channel temperature verification kit (TVK).		
temperature readings.		Call the field service representative to perform a temperature verification test.		
		Perform a temperature verification test using the external multi-channel temperature verification kit (TVK).		
Printing fails.	Printer incorrectly configured on instrument.	Make sure the printer IP address on the instrument matches the IP address from the system administrator.		
	Printer not connected to network.	Connect Ethernet cable to printer.		
	Instrument not connected to	Connect Ethernet cable to instrument.		
	network.	Confirm network is functioning.		
	Ethernet cable bad.	Try a different Ethernet cable.		
		Confirm network is functioning.		
When running on multiple	Instrument(s) not connected to	Connect Ethernet cables to the instruments.		
instruments, can't see instruments in Select	network.	Connect Ethernet cable to the network switch.		
Instruments screen	Switch powered off.	Switch power on.		
	Ethernet cable bad.	Try a different Ethernet cable.		
	IP address conflict.	Change the IP addresses on the instruments.		
	Allow Remote Runs setting in Instrument Name screen set to No.	In Instrument Name screen, change Allow Remote Runs to Yes.		
No screen display	Blown fuse.	Check fuses.		
Screen display missing characters or is illegible.	LCD screen failure.	Request service.		
No response when you power	Not connected to power source.	Switch power on.		
on the instrument.		Connect power cord.		

Observation	Possible cause	Recommended action		
Sample tubes melt during run.	Incompatible consumables used.	Use recommended consumables.		
	Heated cover or sample block(s) too hot.	Run the Heated Cover Verification Test.		
Heating cover is stuck.	Cam, platen, or cover assembly malfunctioning	Request service.		
Sample tubes are crushed or deformed after run.	Incompatible consumables used, or proper trays/retainer accessories not used.	Use recommended consumables.		
Power failure during a run	Power failure during a run	If a power failure occurs during a run, the instrument attempts to continue the run when the power returns. The run begins at the first step of the stage where the power failed. If the time between the initial power failure and the time the power returns is more than 12 hours, the run is not continued.		
		A message stating <i>This run was interrupted</i> by power failure on [date/time] and recovered on [date/time] is displayed upon such an occurance.		
	User did not press Done after completing a run before turning the power off on the instrument.	Always press Done after completing a PCR run on the instrument.		

Return an instrument for service

Before returning an instrument for service, back up the run methods and instrument settings, as described in "Manage Folders" on page 33 and "Recommended instrument settings" on page 53. This practice ensures that you can restore the run methods and other settings in the event that the service center replaces a component that restores the default factory settings, or sends you a replacement instrument.

To return an instrument for service:

- **1.** Contact your local customer care center or technical support group to obtain a copy of the Certificate of Instrument Decontamination, a service notification, a service call number, and, if required, packaging.
- 2. Decontaminate the instrument.

IMPORTANT! If the sample block(s) become contaminated with radioactivity, use a commercially available decontaminant to remove the contamination. If the block(s) cannot be decontaminated, the instrument cannot be returned for service.

- **3.** Complete and sign a copy of the Certificate of Instrument Decontamination.
- **4.** Attach a copy of the Certificate of Instrument Decontamination to the outside of the box, place another copy inside the box, and email a copy to the customer care center.
- **5.** Pack the instrument in the provided packaging, without any accessories or power cords. Include a hard copy of the Certificate of Instrument Decontamination in the box.

Note: DO NOT pack the instrument in packaging not designed for thermal cycler shipment as damage may incur. Repairs for instruments without the Certificate of Instrument Decontamination will be delayed.

6. Affix the provided shipping label to the box, then ship to the designated facility. The repair process requires 1 to 3 weeks, depending on regional location.



Recommended instrument settings

About Instrument

Select **Settings** • **About Instrument** to find out more information about the instrument (e.g., firmware version and instrument statistics).

- Select EULA to view the End User License Agreement. The EULA can be downloaded to a USB drive.
- Select Check updates to find out if updates are available for the instrument .



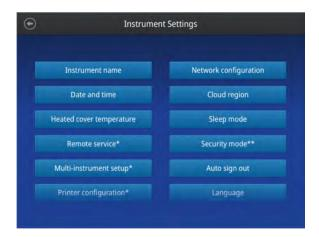
Recommended instrument settings

Select ② (Settings) and access the Settings screen to configure the instrument.



Instrument settings

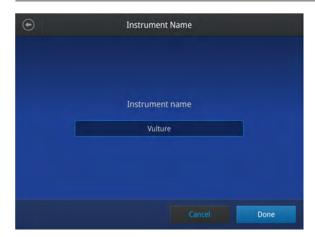
Select Instrument Settings to set the following instrument parameters.



Instrument name

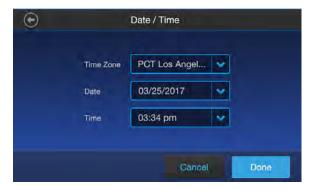
Select the **Instrument name** field to activate the text editor. Enter up to 25 alphanumeric characters to identify the instrument.

Note: The instrument name cannot have spaces. Separate consecutive characters with a hyphen or underscore; for example, *My_Instrument*.



Date and time

- Select the **Time Zone** field to set the time zone.
- Select the **Date/Format** field to choose the date format and set the date.
- Select the Time/Format field to activate the numeric editor to set the time.



· Heated cover temperature

Use the **Off** and **On** toggle to disable or enable activation of the heated cover in idle mode. In the 'On' mode, press the **Edit Temperature** field to activate the numeric editor to set the cover temperature.

Note: The heated cover temperature must be between 30°C and 110°C.



Remote service

Use the **Off** and **On** toggle to disable or enable periodic upload of instrument data to **Connect**.

Multi-instrument setup

Use the **Off** and **On** toggles to make the instrument hidden or visible to other instruments on the network, and to ignore or find other instruments on the network.





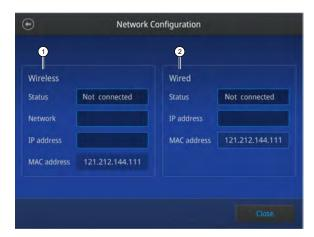
Printer configuration

Select the **Remote Printer IP Address** field to activate the numeric editor, then enter the IP address of the printer to be connected to the instrument.



Network configuration

Select the type of network connection that will be used to connect the instrument to the Internet. For details on using the Wireless and Ethernet options, see "Set up the wired connection" on page 14 and "Set up the wireless connection" on page 17.



- 1 Wireless panel
- (2) Wired panel

• Cloud region (Administrator profile only)

Select the appropriate field to set the cloud region for the instrument.

Sleep mode

Use the **Off** and **On** toggle to disable or enable sleep mode. In the 'On' mode, select the **Edit Time** field to activate the numeric editor to set the time after which the instrument will go from idle mode to standby mode.



• Security mode (Administrator profile only)

Use the toggle to set user accessibility for the instrument.

- Select **Guest accessible** to enable guest access for most instrument functions.
- Select **Login required** to restrict use to those with user accounts.

• Auto sign out (Administrator profile only)

Use the toggle to enable/disable automatic sign out of a user when no activity is detected on the instrument for a selected period of time.

Brightness settings

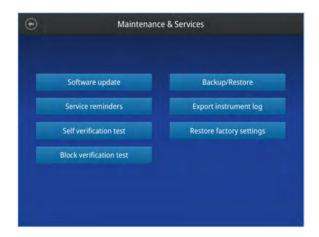
Use the slider to adjust screen brightness.

Language

- Select Settings ➤ Instrument settings ➤ Language ➤ English to enable English language user interface
- Select **Settings ▶ Instrument settings ▶ Language ▶** 简体中文 to enable Chinese language user interface.

Maintenance and services

Select Maintenance & Services to set the following instrument parameters.



- Select **Software Update** to update the System firmware. See "Upgrade the system firmware" on page 46 for instructions on updating the firmware.
- Select **Service reminders** to enable reminders for service after a selected interval of time. Set the time interval by selecting the **Service reminded provided every** field to set the time interval.
- Select Self Verification Test to conduct a check on the instrument hardware. The check includes
 testing the block, heated cover, and other components. See Chapter 5, "Maintenance" for
 instructions on conducting the self-verification test.
- Select Block Verification Test to perform a block verification test. Ensure that you have the Temperature Verification Kit (TVK) (Cat. No. 4377669) before performing this test.
 - Select Verify Block Temperature to carry out the block temperature test. Select the test type, connect the TVK, and insert the TVK probe into a specific zone depending on the test type.
 - Select Verify Cycle Performance to check the cycle performance of the instrument.
- Select Backup/Restore (Administrator profile only) is useful in event of a hardware failure or while setting up multiple instrument runs. See page 19 for details on logging in using the administrator profile.
 - Select Backup Instrument to back up the instrument settings, user accounts, and methods on the instrument to a USB memory device.
 - Select Restore a Backup to view instrument files that are backed up on a USB memory device.
- Select Export Instrument Log to export the instrument logs to a USB memory device. Insert the USB memory device into the USB port before using this feature.
- Select Restore factory settings is used to reset the instrument to the original factory settings
 See "Restore factory settings (Administrator profile only)" on page 49. See page 19 for details on
 logging in using the administrator profile.

Run history

Select **Run History** to display the entire list of runs performed by the instrument.

• Select a particular Run ID to view the details of that run.



• Select **Export** to save the run details to a USB memory device, or **Print** to print the run details.



Note: If you are signed in with the Administrator profile, you can use the **Manage** function to perform the following functions.

- · Delete a run report
- · Select all run reports
- · Export a run report



USB shortcuts

Press **USB Shortcuts** to manage the shortcuts available when a USB memory device is inserted into the instrument.

- Select the shortcuts to be displayed when a USB memory device is inserted.
- Select Show Shortcuts to display the selected shortcuts when a USB memory device is inserted.



Note: Backup/Restore is only available when signed in as administrator.

When a USB memory device is plugged into the USB port, a pop-up dialog box with the pre-selected options is displayed.





Instrument specification and layout

Site requirements

The instrument is for indoor use. Ensure that the installation site:

- Meets the spatial and weight requirements (see "Specifications")
- Meets environmental requirements (see "Environmental requirements")
- Is within 1 m (3 ft.) of an AC power source receptacle
- Is away from water

Environmental requirements

Temperature and humidity requirements

Ensure that the installation site is maintained under the following conditions:

Table 4 Temperature and humidity requirements

Condition	Acceptable range		
Temperature (acceptable range)	15–30°C (59 to 86°F)		
Humidity (acceptable range)	15-80% Relative Humidity, non-condensing		

Avoid placing the instrument adjacent to heaters, cooling ducts, or in direct sunlight. Fluctuations between day and night temperatures can cause system instability. Place away from any equipment that vibrates, such as a refrigerator or centrifuge.

Pollution

The instrument has a Pollution Degree rating of 2. It may be installed in an environment that has non-conductive pollutants only, such as dust particles or wood chips. Typical environments with a Pollution Degree II rating are laboratories and sales and commercial areas.

Altitude

The safety of instrument use was tested for altitudes up to 6000 ft.

Specifications

Technical specifications

Factoria	Specification				
Feature	96-well block	384-well block			
Maximum block ramp rate [1]	6.0°C/sec	5.0°C/sec			
Maximum sample ramp rate	4.4°C/sec	3.5°C/sec			
Temperature accuracy	±0.25°C (35–99.9°C)	±0.25°C (35–99.9°C)			
Temperature range for protocol run	0.0-100.0°C	0.0-100.0°C			
Temperature non-uniformity [2]	≤0.50	≤0.50			
VeriFlex [™] block	 6 VeriFlex[™] Zones Supported: 25°C (5°C zoneto-zone) Allowable ^[3]: 30°C (10°C zone-to-zone) 	_			
PCR volume range	 Supported: 10–100 μL Allowable: 1–100 μL 	 Supported: 5–20 μL Allowable: 1–30 μL 			
Instrument memory	USB, 16GB on-board	USB, 16GB on-board			
Display interface	8 inch color TFT LCD	8 inch color TFT LCD			
IMPORTANT! If the supplied power fluctuates ±10% beyond the rated voltage, a power line regulator may be required. High or low voltages can adversely affect the electronic components of this instrument.	100–120 V ±10%, 50/60 Hz; 200–240 V ±10%, 50/60 Hz	100–120 V ±10%, 50/60 Hz; 200–240 V ±10%, 50/60 Hz			
Installation category	Overvoltage Cat II (Portable equipment)	Overvoltage Cat II (Portable equipment)			
Power consumption	Maximum consumption: 700 W	Maximum consumption: 700 W			
	Average during thermal cycling: 260 W	Average during thermal cycling: 260 W			
	During sleep/ standby: 25 W	During sleep/ standby: 25 W			

 $^{^{[1]}}$ Reaction volume of 1 μ L.

^{[2] 30} seconds after clock starts.

 $^{^{[3]}}$ Temperature accuracy ± 0.50 and Temperature non-uniformity $\leq\!0.75.$

Physical dimensions

Parameter	Instrument footprint	Recommended clearance	Total
Height	21.7 cm (8.54 in)	24 cm (9.44 in) ^[1]	46 cm (18.11 in)
Width	24.5 cm (9.65 in)	0 cm (0 in)	24.5 cm (9.65 in)
Depth	46.5 cm (18.31 in)	23.5 cm (9.25 in) ^[2]	70 cm (27.56 in)
Weight (96-well block model)	12 kg (26.5 lb)		
Weight (384-well block model)	13 kg (28.7 lb)		

^[1] With the heated cover open.

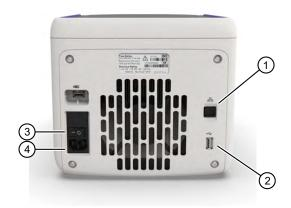
Power and communication port symbols

Port	Description
~	AC power cable port
뀸	10/100 Fast Ethernet port for connecting to a network
•	USB v2.0 port for connecting to an external network drive, jump drive, or other USB storage device
H	USB v2.0 port for connecting to USB-enabled Wi-Fi Card

 $^{^{[2]}}$ Allow enough ventilation space at the rear of the instrument to ensure adequate airflow and cooling.

Location of power point and ports on the instrument

The following graphic displays the location of the power cord insertion point as well as the various ports provided in the instrument.



- 1 Ethernet port
- 2 USB port
- (3) Main power switch
- 4 Power cable port



Predefined Methods

Predefined run methods

The following default methods (distributed across ten categories) are pre-programmed on the instrument. A particular run method can appear in more than one category. These run methods are based on our recommended protocols and should be considered as starting points for the different applications. Your results may vary between instrument models, so some optimization may be needed.

The following table lists the categories and the run methods they include.

Table 5 Pre-defined categories and run methods

Category	Method name
Blank Template	General PCR Run
	Blank PCR 1
	Blank PCR 2
Basic PCR	• AmpliTaq [™] 360
	 AmpliTaq[™]
	Basic PCR Taq
	 DreamTaq[™] DNA Polymerase 3-Step
Hot Start	AmpliTaq Gold [™] 360
	 AmpliTaq Gold[™]
	 AmpliTaq Gold[™] Fast PCR Run
	 DreamTaq[™] Hot Start DNA Master Mix
	 Phusion[™] Hot Start II PCR Master Mix 2-Step
	 Phusion[™] Hot Start II PCR Master Mix 3-Step
	 Platinum[™] Direct PCR Universal Master Mix
	 Platinum[™] Taq DNA Polymerase
	Platinum [™] Taq Master Mix
	 Platinum[™] SuperFi II PCR Master Mix
	 Platinum[™] SuperFi II PCR Master Mix 2-Step
	 Platinum[™] SuperFi II PCR Master Mix 3-Step
	 Platinum[™] II Taq Hot-Start PCR Master Mix 2-Step
	 Platinum[™] II Taq Hot-Start PCR Master Mix 3-Step

Table 5 Pre-defined categories and run methods (continued)

Category	Method name
Sequencing	 BigDye[™] Direct Cycle Sequence BigDye[™] Direct PCR Run BigDye[™] Kit Fast Run BigDye[™] Kit Standard Run Collibri[™] ES DNA Prep Collibri[™] PS DNA Prep DreamTaq[™] Hot Start DNA Master Mix Phusion[™] Hot Start II PCR Master Mix 2-Step Phusion[™] Hot Start II PCR Master Mix 3-Step Platinum[™] Direct PCR Universal Master Mix Platinum[™] SuperFi PCR Master Mix 2-Step Platinum[™] SuperFi PCR Master Mix 3-Step Platinum[™] SuperFi II PCR Master Mix Platinum[™] Taq Master Mix Platinum[™] Taq Master Mix Platinum[™] II Taq Hot-Start PCR Master Mix 2-Step Platinum[™] II Taq Hot-Start PCR Master Mix 3-Step
Optimization	 Time Release Run Touchdown PCR Run DreamTaq[™] Hot Start DNA Master Mix Phusion[™] Hot Start II PCR Master Mix 2-Step Phusion[™] Hot Start II PCR Master Mix 3-Step Platinum[™] Direct PCR Universal Master Mix Platinum[™] SuperFi PCR Master Mix 2-Step Platinum[™] SuperFi PCR Master Mix 3-Step Platinum[™] SuperFi II PCR Master Mix Platinum[™] Taq Master Mix Platinum[™] Taq Hot-Start PCR Master Mix 2-Step Platinum[™] II Taq Hot-Start PCR Master Mix 3-Step
RT-PCR	SuperScript [™] III 1-Step System SuperScript [™] IV One-Step RT-PCR System

Table 5 Pre-defined categories and run methods (continued)

Category	Method name
High Fidelity	 AccuPrime[™] PFX 2-Step AccuPrime[™] PFX 3-Step AccuPrime[™] Taq High Fidelity Phusion[™] Flash PCR Master Mix 2-Step Phusion[™] Flash PCR Master Mix 3-Step Phusion[™] High Fidelity DNA Polymerase 2-Step Phusion[™] High Fidelity DNA Polymerase 3-Step Phusion[™] Hot Start II PCR Master Mix 2-Step Phusion[™] Hot Start II PCR Master Mix 3-Step Platinum[™] Taq High Fidelity Platinum[™] SuperFi II PCR Master Mix Platinum[™] SuperFi II PCR Master Mix 2-Step Platinum[™] SuperFi II PCR Master Mix 2-Step Platinum[™] SuperFi II PCR Master Mix 3-Step
High Specificity	 AccuPrime[™] PFX 2-Step AccuPrime[™] Taq DNA Polymerase AccuPrime[™] Taq High Fidelity DreamTaq[™] Hot Start DNA Master Mix Phusion[™] Hot Start II PCR Master Mix 2-Step Phusion[™] Hot Start II PCR Master Mix 3-Step Platinum[™] Direct PCR Universal Master Mix Platinum[™] SuperFi PCR Master Mix 2-Step Platinum[™] SuperFi PCR Master Mix 3-Step Platinum[™] SuperFi II PCR Master Mix Platinum[™] Taq Master Mix Platinum[™] Taq Master Mix Platinum[™] II Taq Hot-Start PCR Master Mix 3-Step Platinum[™] II Taq Hot-Start PCR Master Mix 3-Step
Next-Gen Sequencing	TargetSeq [™] Exome Enrichment
Long PCR	 AccuPrime[™] Taq High Fidelity Phusion[™] Hot Start II PCR Master Mix 3-Step Platinum[™] SuperFi PCR Master Mix Platinum[™] SuperFi II PCR Master Mix Platinum[™] Taq High Fidelity

These run methods cannot be edited or deleted, but if you save them with a different name or to a different folder, you can edit them.

The following tables describe the method details.

Note: For illustration purpose, the tables display maximum ramp rate.

AccuPrime[™] PFX 2-Step Method

Cycle/Step		Stage				Reaction	Cover
		1	2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	94.0	94.0	4.0			
	Hold Time	120	15	Infinite			
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		68.0				
	Hold Time		60				

 $^{^{[1]}~~\}text{Values for VeritiPro}^{^{\text{\tiny TM}}}~\text{Thermal Cycler}.$

AccuPrime[™] PFX 3-Step Method

Cycle/Step		Stage				Reaction	Cover
		1	2	3	4	volume (µL)	temp. (°C)
Cycles		1	25	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	94.0	94.0	4.0			
	Hold Time	120	15	Infinite			
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		55.0				
	Hold Time		15				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		68.0				
	Hold Time		60				

 $^{^{[1]}~~\}text{Values for VeritiPro}^{^{\intercal}}~\text{Thermal Cycler}.$

AccuPrime[™] Taq DNA Polymerase Method

Cycle/Step			Sta	age		Reaction volume (µL)	Cover temp. (°C)
		1	2	3	4		
Cycles		1	25	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	94.0	94.0	4.0			
	Hold Time	120	15	Infinite			
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		55.0				
	Hold Time		15				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		68.0				
	Hold Time		60				

^[1] Values for VeritiPro[™] Thermal Cycler.

AccuPrime[™] Taq High Fidelity Method

Cycle/Step			Sta	Reaction	Cover		
		1	2	3	4	volume (µL)	temp. (°C)
Cycles		1	25	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	94.0	94.0	4.0			
	Hold Time	15	15	Infinite			
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		52.0				
	Hold Time		15				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		68.0				
	Hold Time		60				

^[1] Values for VeritiPro[™] Thermal Cycler.

AmpliTaq[™] 360 Method

Ovela/Otem			Sta	age		Reaction Cov			
	Cycle/Step		2	3	4	volume (µL)	temp. (°C)		
Cycles		1	35	1	_	50	105.0		
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_				
	Temperature	94.0	95.0	72.0					
	Hold Time	180	30	420					
Step 2	Ramp Rate	_	6.0	6.0	_				
	Temperature		58.0	4.0					
	Hold Time		30	Infinite					
Step 3	Ramp Rate	_	6.0	_	_				
	Temperature		72.0						
	Hold Time		60						

^[1] Values for VeritiPro[™] Thermal Cycler.

AmpliTaq Gold[™] 360 Method

Cycle/Step			Sta	age		Cover	
		1	2	3	4	volume (μL)	temp. (°C)
Cycles		1	35	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	95.0	95.0	72.0			
	Hold Time	300	15	420			
Step 2	Ramp Rate	_	6.0	6.0	_		
	Temperature		55.0	4.0			
	Hold Time		30	Infinite			
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		60				

^[1] Values for VeritiPro[™] Thermal Cycler.

AmpliTaq Gold[™] Fast PCR Method

Cycle/Step			Sta	age		Reaction	Cover
		1	2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	95.0	96.0	72.0			
	Hold Time	600	3	10			
Step 2	Ramp Rate	_	6.0	6.0	_		
	Temperature		62.0	4.0			
	Hold Time		3	Infinite			
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		68.0				
	Hold Time	1	5]			

^[1] Values for VeritiPro[™] Thermal Cycler.

AmpliTaq Gold[™] Method

Cycle/Step			Sta	Reaction	Cover		
		1	2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	95.0	95.0	72.0			
	Hold Time	300	15	420			
Step 2	Ramp Rate	_	6.0	6.0	_		
	Temperature		55.0	4.0			
	Hold Time		15	Infinite			
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		30				

^[1] Values for VeritiPro[™] Thermal Cycler.

AmpliTaq[™] Method

Cycle/Step			Sta	age		Reaction	Cover temp. (°C)
		1	2	3	4	volume (µL)	
Cycles		1	35	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	95.0	95.0	72.0			
	Hold Time	60	15	420			
Step 2	Ramp Rate	_	6.0	6.0	_		
	Temperature		55.0	4.0			
	Hold Time		15	Infinite			
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		30				

^[1] Values for VeritiPro[™] Thermal Cycler.

Basic PCR Taq Method

Cycle/Step			Sta	age		Cover	
		1	2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	95.0	95.0	72.0			
	Hold Time	60	15	420			
Step 2	Ramp Rate	_	6.0	6.0	_		
	Temperature		55.0	4.0			
	Hold Time		15	Infinite			
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		30				

^[1] Values for VeritiPro[™] Thermal Cycler.

BigDye[™] Direct Cycle Sequence Method

0	Cycle/Step			Stage			Reaction	Cover
Су			2	3	4	5	volume (µL)	temp. (°C)
Cycles		1	1	1	25	1	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	6.0	6.0		
	Temperature	37.0	80.0	96.0	96.0	4.0		
	Hold Time	900	120	60	10	Infinite		
Step 2	Ramp Rate	_	_	_	6.0	_		
	Temperature				50.0			
	Hold Time				5			
Step 3	Ramp Rate	_	_	_	6.0	_		
	Temperature				60.0			
	Hold Time				75			

^[1] Values for VeritiPro[™] Thermal Cycler.

BigDye[™] Direct PCR Run Method

			Sta	Reaction	Cover		
	Cycle/Step		2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	95.0	96.0	72.0			
	Hold Time	600	3	120			
Step 2	Ramp Rate	_	6.0	6.0	_		
	Temperature		62.0	4.0			
	Hold Time		15	Infinite			
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		68.0				
	Hold Time		30				

 $^{^{[1]}}$ Values for VeritiPro $^{™}$ Thermal Cycler.

BigDye[™] Kit Fast Method

			Sta	Reaction	Cover		
	Cycle/Step		2	3	4	volume (µL)	temp. (°C)
Cycles		1	25	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	96.0	96.0	4.0			
	Hold Time	60	10	Infinite			
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		50.0				
	Hold Time		5				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		60.0				
	Hold Time		75				

^[1] Values for VeritiPro[™] Thermal Cycler.

BigDye[™] Kit Standard Method

0	vola/Stan		Sta	Reaction	Cover		
Cycle/Step		1	2	3	4	volume (µL)	temp. (°C)
Cycles		1	25	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	96.0	96.0	4.0			
	Hold Time	60	10	Infinite			
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		50.0				
	Hold Time		5				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		60.0				
	Hold Time		240				

^[1] Values for VeritiPro[™] Thermal Cycler.

Collibri[™] ES DNA Prep Method

Ovele/Stern			Sta	Reaction	Cover		
	Cycle/Step		2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	94.0	94.0	4.0			
	Hold Time	120	15	Infinite			
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		68.0				
	Hold Time		60				

 $^{^{[1]}~}$ Values for VeritiPro $^{^{\text{\tiny{TM}}}}$ Thermal Cycler.

Collibri[™] PS DNA Prep Method

	l.a /Ota		Sta	Reaction	Cover		
	Cycle/Step		2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	94.0	94.0	4.0			
	Hold Time	120	15	Infinite			
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		55.0				
	Hold Time		30				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		68.0				
	Hold Time		60				

 $^{^{[1]}~~\}text{Values for VeritiPro}^{^{\text{\tiny{TM}}}}~\text{Thermal Cycler}.$

DreamTaq[™] Hot Start PCR Master Mix Method

0.	Cycle/Step		Sta	Reaction	Cover		
C			2	3	4	volume (µL)	temp. (°C)
Cycles		1	25–40	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	95.0	95.0	72.0			
	Hold Time	120	30	600			
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		60.0				
	Hold Time		60				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		60				

^[1] Values for VeritiPro[™] Thermal Cycler.

DreamTaq[™] DNA Polymerase 3-Step

	Cycle/Step		Sta	Reaction	Cover		
Сусіе/Зіер		1	2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	1	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	6.0		
	Temperature	95.0	95.0	72.0	4.0		
	Hold Time	120	30	600	Infinite		
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		60.0				
	Hold Time		30				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		60				

 $^{^{[1]}}$ Values for VeritiPro $^{\text{\tiny{TM}}}$ Thermal Cycler.

General PCR Method

			Sta	Reaction	Cover		
	Cycle/Step		2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	95.0	95.0	72.0			
	Hold Time	60	15	420			
Step 2	Ramp Rate	_	6.0	6.0	_		
	Temperature		55.0	4.0			
	Hold Time		15	Infinite			
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time	1	30				

^[1] Values for VeritiPro[™] Thermal Cycler.

Phusion[™] Flash PCR Master Mix 2-Step

0	vala/Stan		Sta	Reaction	Cover		
C	Cycle/Step		2	3	4	volume (µL)	temp. (°C)
Cycles		1	30	1	1	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	6.0		
	Temperature	98.0	98.0	72.0	4.0		
	Hold Time	10	1	60	Infinite		
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		15				
Step 3	Ramp Rate	_	_	_	_		
	Temperature						
	Hold Time						

^[1] Values for VeritiPro[™] Thermal Cycler.

Phusion[™] Flash PCR Master Mix 3-Step

			Sta	Reaction	Cover		
	Cycle/Step		2	3	4	volume (µL)	temp. (°C)
Cycles		1	30	1	1	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	6.0		
	Temperature	98.0	98.0	72.0	4.0		
	Hold Time	10	1	60	Infinite		
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		65.0				
	Hold Time		5				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		15				

^[1] Values for VeritiPro[™] Thermal Cycler.

Phusion[™] High Fidelity DNA Polymerase 2-Step

0	Cycle/Step		Sta	Reaction	Cover		
C			2	3	4	volume (µL)	temp. (°C)
Cycles		1	30	1	1	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	6.0		
	Temperature	98.0	98.0	72.0	4.0		
	Hold Time	30	10	600	Infinite		
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		30				
Step 3	Ramp Rate	_	_	_	_		
	Temperature						
	Hold Time						

^[1] Values for VeritiPro[™] Thermal Cycler.

Phusion[™] High Fidelity DNA Polymerase 3-Step

			Sta	Reaction	Cover		
G	ycle/Step	1	2	3	4	volume (µL)	temp. (°C)
Cycles		1	30	1	1	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	6.0		
	Temperature	98.0	98.0	72.0	4.0		
	Hold Time	30	10	600	Infinite		
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		65.0				
	Hold Time		30				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		30				

^[1] Values for VeritiPro[™] Thermal Cycler.

Phusion[™] Hot Start II PCR Master Mix 2-Step

	Cycle/Step		Sta	Reaction	Cover		
C			2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	1	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	6.0		
	Temperature	98.0	98.0	72.0	4.0		
	Hold Time	30	10	300	Infinite		
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		30				
Step 3	Ramp Rate	_	_	_	_		
	Temperature						
	Hold Time						

^[1] Values for VeritiPro[™] Thermal Cycler.

Phusion[™] Hot Start II PCR Master Mix 3-Step

0.			Sta	Reaction	Cover		
C	Cycle/Step		2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	1	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	6.0		
	Temperature	98.0	98.0	72.0	4.0		
	Hold Time	30	10	300	Infinite		
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		60.0				
	Hold Time		10				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		30				

^[1] Values for VeritiPro[™] Thermal Cycler.

Platinum[™] Direct PCR Universal Master Mix Method

			Sta	Reaction	Cover		
C	Cycle/Step		2	3	4	volume (µL)	temp. (°C)
Cycles		1	40	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	94.0	94.0	4.0			
	Hold Time	120	15	Infinite			
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		60.0				
	Hold Time		15				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		68.0				
	Hold Time		20				

^[1] Values for VeritiPro[™] Thermal Cycler.

Platinum[™] SuperFi PCR Master Mix Method

			Sta	Reaction	Cover		
	Cycle/Step		2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	1	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	6.0		
	Temperature	95.0	95.0	68.0	4.0		
	Hold Time	120	10	300	Infinite		
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		60.0				
	Hold Time		10				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		68.0				
	Hold Time	1	30				

^[1] Values for VeritiPro[™] Thermal Cycler.

Platinum[™] SuperFi PCR Master Mix 2-Step

0	vala/Stan		Sta	Reaction	Cover		
C	Cycle/Step		2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	1	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	6.0		
	Temperature	98.0	98.0	72.0	4.0		
	Hold Time	30	10	300	Infinite		
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		30				
Step 3	Ramp Rate	_	_	_	_		
	Temperature						
	Hold Time						

^[1] Values for VeritiPro[™] Thermal Cycler.

Platinum[™] SuperFi PCR Master Mix 3-Step

0.			Sta	Reaction	Cover		
C	Cycle/Step		2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	1	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	6.0		
	Temperature	98.0	98.0	72.0	4.0		
	Hold Time	30	10	300	Infinite		
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		60.0				
	Hold Time		10				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		30				

^[1] Values for VeritiPro[™] Thermal Cycler.

Platinum[™] SuperFi II PCR Master Mix Method

0	rala/Stan		Sta	Reaction	Cover		
C ₁	Cycle/Step		2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	1	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	6.0		
	Temperature	98.0	98.0	72.0	4.0		
	Hold Time	30	10	300	Infinite		
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		60.0				
	Hold Time		10				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		30				

^[1] Values for VeritiPro[™] Thermal Cycler.

Platinum[™] Taq DNA Polymerase Method

	Cycle/Step		Sta	Reaction	Cover		
			2	3	4	volume (µL)	temp. (°C)
Cycles		1	30	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	94.0	94.0	4.0			
	Hold Time	30	30	Infinite			
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		55.0				
	Hold Time		30				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		60				

^[1] Values for VeritiPro[™] Thermal Cycler.

Platinum[™] Taq High Fidelity Method

			Sta	Reaction	Cover		
	Cycle/Step		2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	94.0	94.0	4.0			
	Hold Time	30	15	Infinite			
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		55.0				
	Hold Time		15				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		68.0				
	Hold Time		60				

^[1] Values for VeritiPro[™] Thermal Cycler.

Platinum[™] Taq Master Mix Method

	hada (Otan		Sta	Reaction	Cover		
	ycle/Step	1	2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	94.0	94.0	4.0			
	Hold Time	120	30	Infinite			
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		55.0				
	Hold Time		30				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		60				

^[1] Values for VeritiPro[™] Thermal Cycler.

Platinum[™] II Taq Hot-Start PCR Master Mix 2-Step

	David & 101 and		Sta	Reaction	Cover		
Cycle/Step		1	2	3	4	volume (µL)	temp. (°C)
Cycles		1	30	1	1	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	6.0		
	Temperature	98.0	98.0	72.0	4.0		
	Hold Time	30	10	600	Infinite		
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		30				
Step 3	Ramp Rate	_	_	_	_		
	Temperature						
	Hold Time						

 $^{^{[1]}}$ Values for VeritiPro $^{^{\text{\tiny TM}}}$ Thermal Cycler.

Platinum[™] II Taq Hot-Start PCR Master Mix 3-Step

	Cycle/Step		Sta	age	Reaction	Cover	
			2	3	4	volume (µL)	temp. (°C)
Cycles		1	35	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	94.0	94.0	4.0			
	Hold Time	120	15	Infinite			
Step 2	Ramp Rate	_	6.0	_	_		
	Temperature		60.0				
	Hold Time		15				
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		68.0				
	Hold Time		15				

^[1] Values for VeritiPro[™] Thermal Cycler.

SuperScript[™] III 1-Step System Method

0	Cycle/Step		Stage				Cover
C			2	3	4	volume (µL)	temp. (°C)
Cycles		1	40	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	60.0	94.0	68.0			
	Hold Time	900	15	300			
Step 2	Ramp Rate	6.0	6.0	6.0	_		
	Temperature	94.0	55.0	4.0			
	Hold Time	120	30	Infinite			
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		68.0				
	Hold Time		60				

^[1] Values for VeritiPro[™] Thermal Cycler.

SuperScript[™] IV One-Step RT-PCR System Method

	Cycle/Step		Sta	age		Reaction	Cover
			2	3	4	volume (µL)	temp. (°C)
Cycles		1	40	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	60.0	94.0	68.0			
	Hold Time	900	15	300			
Step 2	Ramp Rate	6.0	6.0	6.0	_		
	Temperature	94.0	55.0	4.0			
	Hold Time	120	30	Infinite			
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		68.0				
	Hold Time		60				

^[1] Values for VeritiPro[™] Thermal Cycler.

Time Release Method

0-	Cycle/Step		Sta	age		Reaction	Cover
C			2	3	4	volume (µL)	temp. (°C)
Cycles		1	40	1	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0	6.0	_		
	Temperature	95.0	95.0	72.0			
	Hold Time	60	15	420			
Step 2	Ramp Rate	_	6.0	6.0	_		
	Temperature		55.0	4.0			
	Hold Time		15	Infinite			
Step 3	Ramp Rate	_	6.0	_	_		
	Temperature		72.0				
	Hold Time		30				

^[1] Values for VeritiPro[™] Thermal Cycler.

TargetSeq[™] Exome Enrichment Method

Cycle/Step			St	age		Reaction	Cover
		1	2	3	4	volume (µL)	temp. (°C)
Cycles		72	_	_	_	50	105.0
Step 1	Ramp Rate ^[1]	6.0	_	_	_		
	Temperature	47.0					
	Hold Time	3600					

 $^{^{[1]}~}$ Values for VeritiPro $^{^{\text{\tiny TM}}}$ Thermal Cycler.

Touchdown PCR Method

Ovela/Otem			Sta	age	Reaction	Cover	
	Cycle/Step		2	3	4	volume (µL)	temp. (°C)
Cycles		20	10			50	105.0
Step 1	Ramp Rate ^[1]	6.0	6.0				
	Temperature	94.0	94.0				
	Hold Time	15	15				
Step 2	Ramp Rate	6.0	6.0				
	Temperature	65.0	55.0	_	_		
	Hold Time	30	30				
	AutoDelta Temperature	-0.5	_				
	AutoDelta Starting Cycle	1	_				

 $^{^{[1]}~~\}text{Values for VeritiPro}^{^{\text{\tiny TM}}}~\text{Thermal Cycler}.$



About the Thermal Cycler Fleet Control Software

The Applied Biosystems[™] Thermal Cycler Fleet Control Software (Cat. No. A40070) is used to manage users, methods, run reports, and multiple PCR instruments through a single user interface. The software is compatible with all Applied Biosystems[™] Thermal Cyclers. The Thermal Cycler Fleet Control Software can be purchased separately, and is not included with the VeritiPro[™] Thermal Cycler. For more information visit thermofisher.com.

Connect the thermal cycler to the network

- 1. Connect the thermal cycler to your network using an Ethernet cable.
- 2. Go to Settings Menu ➤ Admin Menu ➤ Set IP Address to set up your thermal cycler network configuration.



Ordering information

Instrument part numbers

Contact your regional sales office to order kits, accessories, consumables, and spare parts for the Veriti Pro^{TM} 96-well Thermal Cycler (Cat. No. A48141) and Veriti Pro^{TM} 384-well Thermal Cycler (Cat. No. A48140).

Consumables

Product	Cat. No.
Single Tubes	
MicroAmp [™] Reaction Tube with Cap, 0.2 mL	N8010540
	N8011540
MicroAmp [™] Reaction Tube with Cap, 0.2 mL, assorted colors	N8010840
MicroAmp [™] Reaction Tube with Cap, 0.2 mL, autoclaved	N8010612
MicroAmp [™] Reaction Tube without Cap, 0.2 mL	N8010533
	N8011533
MicroAmp [™] Reaction Tube without Cap, 0.2 mL, assorted colors	N8010833
MicroAmp [™] Optical Tube without Cap, 0.2 mL	N8010933
Reaction Trays	
MicroAmp [™] 96-Well Tray for VeriFlex [™] Block	4379983
MicroAmp [™] 96-Well Tray/Retainer Set	4381850
96-Well Plates	
MicroAmp [™] Optical 96-Well Reaction Plate	N8010560
	4316813
MicroAmp [™] Optical 96-Well Reaction Plate with Barcode	4306737
	4326659
MicroAmp [™] Optical 96-Well Reaction Plate with Barcode & Optical Caps	403012

Appendix F Ordering information Consumables

(continued)

Product	Cat. No.
MicroAmp [™] Optical 96-Well Reaction Plate with Barcode & Optical Adhesive Films	4314320
MicroAmp [™] EnduraPlate [™] Optical 96-Well Clear Reaction Plates with Barcode	4483354
	4483352
MicroAmp [™] EnduraPlate [™] Optical 96-Well Blue Reaction Plates with Barcode	4483343
MicroAmp [™] EnduraPlate [™] Optical 96-Well Green Reaction Plates with Barcode	4483349
MicroAmp [™] EnduraPlate [™] Optical 96-Well Red Reaction Plates with Barcode	4483350
MicroAmp [™] EnduraPlate [™] Optical 96-Well Yellow Reaction Plates with Barcode	4483395
MicroAmp [™] EnduraPlate [™] Optical 96-Well Multicolor Reaction Plates with Barcode	4483355
	4483356
MicroAmp [™] EnduraPlate [™] Optical 96-Well Clear GPLE Reaction Plates with Barcode	4483348
	4483351
384-Well Plates	
MicroAmp [™] Optical 384-Well Reaction Plate with Barcode	4309849
MicroAmp [™] EnduraPlate [™] Optical 384-Well Clear Reaction Plates with Barcode	4483285
MicroAmp [™] EnduraPlate [™] Optical 384-Well Yellow Reaction Plates with Barcode	4483315
MicroAmp [™] EnduraPlate [™] Optical 384-Well Clear GPLE Reaction Plates with Barcode	4483319
MicroAmp [™] EnduraPlate [™] Optical 384-Well Blue Reaction Plates with Barcode	4483320
MicroAmp [™] EnduraPlate [™] Optical 384-Well Green Reaction Plates with Barcode	4483321
MicroAmp [™] EnduraPlate [™] Optical 384-Well Red Reaction Plates with Barcode	4483322
8-Well Strips	
MicroAmp [™] 8-Tube Strip, 0.2 mL	N8010580
MicroAmp [™] 8-Tube Strip, 0.2 mL, assorted colors	N8010838
MicroAmp [™] Optical 8-Tube Strip, 0.2 mL	4316567
MicroAmp [™] 8-Cap Strip, clear	N8010535
	N8011535
MicroAmp [™] 8-Cap Strip, assorted colors	N8010835
MicroAmp [™] Optical 8-Cap Strips	4323032
Accessories	
MicroAmp [™] Adhesive Film Applicator	4333183

(continued)

Product	Cat. No.
MicroAmp [™] Clear Adhesive Film	4306311
MicroAmp [™] Multi Removal Tool	4313950
MicroAmp [™] Cap Installing Tool	4330015
MicroAmp [™] 96-Well Base	N8010531
MicroAmp [™] Splash-Free 96-Well Base	4312063
MicroAmp [™] Centrifuge Adapter	N8013822

Safety





WARNING! GENERAL SAFETY. Using this product in a manner not specified in the user documentation may result in personal injury or damage to the instrument or device. Ensure that anyone using this product has received instructions in general safety practices for laboratories and the safety information provided in this document.

- Before using an instrument or device, read and understand the safety information provided in the user documentation provided by the manufacturer of the instrument or device.
- Before handling chemicals, read and understand all applicable Safety Data Sheets (SDSs) and use appropriate personal protective equipment (gloves, gowns, eye protection, and so on). To obtain SDSs, see the "Documentation and Support" section in this document.

Symbols on the instrument

Symbols may be found on the instrument to warn against potential hazards or convey important safety information. In this document, the hazard symbol is used along with one of the following user attention words:

- **CAUTION!** Indicates a potentially hazardous situation that, if not avoided, may result in minor or moderate injury. It may also be used to alert against unsafe practices.
- WARNING! Indicates a potentially hazardous situation that, if not avoided, could result in death or serious injury.
- DANGER! Indicates an imminently hazardous situation that, if not avoided, will result in death or serious injury.

Symbol	English	Français
3 5	Environmental protection symbol of the China RoHS directive. The number in the symbol indicates the "Environment-friendly Use Period" of the product in years. The symbol is used if a substance restricted in China is used in excess of the maximum permitted limit.	Symbole de protection de l'environnement de la directive RoHS de la Chine. Le nombre indiqué dans le symbole indique "période d'utilisation favorable à l'environnement" du produit au cours des années. Le symbole est utilisé si une substance restreinte en Chine est utilisée au-delà de la limite maximale autorisée.

Symbol	English	Français
<u> </u>	Caution, risk of danger Consult the manual for further safety information.	Attention, risque de danger Consulter le manuel pour d'autres renseignements de sécurité.

(continued)

Symbol	English	Français		
	Protective conductor terminal (main ground)	Borne de conducteur de protection (mise à la terre principale)		
	Do not dispose of this product in unsorted municipal waste	Ne pas éliminer ce produit avec les déchets usuels non soumis au tri sélectif.		
	environmental impact from disposal of electronic waste, do not dispose of electronic waste in unsorted municipal waste. Follow local municipal waste ordinances for proper disposal provision and contact customer service for information about responsible disposal options.	mise en garde! Pour minimiser les conséquences négatives sur l'environnement à la suite de l'élimination de déchets électroniques, ne pas éliminer ce déchet électronique avec les déchets usuels non soumis au tri sélectif. Se conformer aux ordonnances locales sur les déchets municipaux pour les dispositions d'élimination et communiquer avec le service à la clientèle pour des renseignements sur les options d'élimination responsable.		

Safety symbols on this instrument

Symbol and description					
CAUTION!	Hot surface.				
CAUTION!	Potential biohazard.				

Location of safety labels on the instrument



Figure 2 Location of the safety label on the rear panel of the instrument



Figure 3 Location of the safety label on the sample block

China RoHS Compliance

本表格依据 SJ/T11364 的规定编制

The following table is prepared in accordance with the provisions of SJ/T11364.

表 6

产品中有害物质的名称及含量									
China EEP Hazardous Substances Information									
部件名称 Component name	VeritiPro [™] 96-well Thermal Cycler (Cat. No. A48141) and VeritiPro [™] 384-well Thermal Cycl (Cat. No. A48140)								
	有害物质								
	Hazardous Substances								
	铅	汞	镉	六价铬	多溴联苯	多溴二苯醚			
	(Pb)	(Hg)	(Cd)	(Cr (VI))	(PBB)	(PBDE)			
电路板	×	0	0	0	0	0			
PCBA's	^	O							
电源供应器	×	0	0	0	0	0			
Power supply		0)			0			
电机组件									
Electromechanical assemblies	×	0	0	0	0	0			

〇: 表示该有害物质在该部件所有均质材料中的含量均在 GB/T26572 规定的限量要求以下.

Indicates that the concentration of the hazardous substance in all homogenous materials for the part is below the relevant threshold of the **GB/T 26572** standard.

×: 表示该有害物质至少在该部件的某一均质材料中的含量超出 GB/T26572 规定的限量要.

Indicates that the concentration of the hazardous substance in at least one homogenous material of the part is above the relevant threshold of **GB/T 26572** standard.

Safety information for instruments not manufactured by Thermo Fisher Scientific

Some of the accessories provided as part of the instrument system are not designed or built by Thermo Fisher Scientific. Consult the manufacturer's documentation for the information needed for the safe use of these products.

Instrument safety

General



CAUTION! Do not remove instrument protective covers. If you remove the protective instrument panels or disable interlock devices, you may be exposed to serious hazards including, but not limited to, severe electrical shock, laser exposure, crushing, or chemical exposure.

Physical injury



CAUTION! Moving Parts. Moving parts can crush, pinch and cut. Keep hands clear of moving parts while operating the instrument. Disconnect power before servicing.



WARNING! Sharp Edges. The edges of this container are sharp.

Electrical safety



WARNING! Ensure appropriate electrical supply. For safe operation of the instrument:

- Plug the system into a properly grounded receptacle with adequate current capacity.
- · Ensure the electrical supply is of suitable voltage.
- Never operate the instrument with the ground disconnected. Grounding continuity is required for safe operation of the instrument.



WARNING! Power Supply Line Cords. Use properly configured and approved line cords for the power supply in your facility.



WARNING! Disconnecting Power. To fully disconnect power either detach or unplug the power cord, positioning the instrument such that the power cord is accessible.



Cleaning and decontamination



CAUTION! Cleaning and Decontamination. Use only the cleaning and decontamination methods specified in the manufacturer's user documentation. It is the responsibility of the operator (or other responsible person) to ensure the following requirements are met:

- No decontamination or cleaning agents are used that could cause a HAZARD as a result of a reaction with parts of the equipment or with material contained in the equipment.
- The instrument is properly decontaminated a) if hazardous material is spilled onto or into the equipment, and/or b) prior to having the instrument serviced at your facility or sending the instrument for repair, maintenance, trade-in, disposal, or termination of a loan (decontamination forms may be requested from customer service).
- Before using any cleaning or decontamination methods (except those recommended by the manufacturer), users should confirm with the manufacturer that the proposed method will not damage the equipment.

Safety and electromagnetic compatibility (EMC) standards

The instrument design and manufacture complies with the following standards and requirements for safety and electromagnetic compatibility.

Safety

Reference	Description	
EU Directive 2014/35/EU	European Union "Low Voltage Directive"	
IEC 61010-1 EN 61010-1 UL 61010-1 CAN/CSA C22.2 No. 61010-1	Safety requirements for electrical equipment for measurement, control, and laboratory use – Part 1: General requirements	
IEC 61010-2-010 EN 61010-2-010 UL 61010-1-2-010 CAN/CSA C22.2 No. 61010-2-010	Safety requirements for electrical equipment for measurement, control and laboratory use – Part 2-010: Particular requirements for laboratory equipment for the heating of materials	
IEC 61010-2-081 EN 61010-2-081 UL 61010-1-2-081 CAN/CSA C22.2 No. 61010-2-081	Safety requirements for electrical equipment for measurement, control and laboratory use – Part 2-081: Particular requirements for automatic and semi-automatic laboratory equipment for analysis and other purposes	

EMC

Reference	Description		
Directive 2014/30/EU	European Union "EMC Directive"		
IEC 61326-1 EN 61326-1	Electrical Equipment for Measurement, Control and Laboratory Use – EMC Requirements – Part 1: General Requirements		
AS/NZS CISPR11	Limits and Methods of Measurement of Electromagnetic Disturbance Characteristics of Industrial, Scientific, and Medical (ISM) Radiofrequency Equipment		
ICES-001, Issue 4	Industrial, Scientific and Medical (ISM) Radio Frequency Generators		
FCC Part 15 Subpart B (47 CFR)	U.S. Standard Radio Frequency Devices		

Class A notice

This equipment has been tested and found to comply with the limits for a Class A digital device, pursuant to part 15 of the FCC Rules. These limits are designed to provide reasonable protection against harmful interference when the equipment is operated in a commercial environment. This equipment generates, uses, and can radiate radio frequency energy and, if not installed and used in accordance with the instruction manual, may cause harmful interference to radio communications. Operation of this equipment in a residential area is likely to cause harmful interference in which case the user will be required to correct the interference at his own expense.

Environmental design standards

Reference	Description
EU Directive 2012/19/EU	European Union "WEEE Directive" - Waste electrical and electronic equipment
EU Directive 2011/65/EU Commission Delegated Directive (EU) 2015/863	European Union "RoHS Directive" — Restriction of hazardous substances in electrical and electronic equipment
GB/T 26572-2011	RRequirements of concentration limits for certain restricted substances in electrical and electronic products
SJ/T 11364-2014	Marking for the restricted use of hazardous substances in electronic and electrical products

Chemical safety



WARNING! GENERAL CHEMICAL HANDLING. To minimize hazards, ensure laboratory personnel read and practice the general safety guidelines for chemical usage, storage, and waste provided below. Consult the relevant SDS for specific precautions and instructions:

- Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. To obtain SDSs, see the "Documentation and Support" section in this document.
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing).
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood).
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer's cleanup procedures as recommended in the SDS.
- · Handle chemical wastes in a fume hood.
- Ensure use of primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary container.
 Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)
- After emptying a waste container, seal it with the cap provided.
- Characterize (by analysis if necessary) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure that the waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.
- **IMPORTANT!** Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.

Appendix G Safety Biological hazard safety

Biological hazard safety



WARNING! BIOHAZARD. Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. Conduct all work in properly equipped facilities with the appropriate safety equipment (for example, physical containment devices). Safety equipment can also include items for personal protection, such as gloves, coats, gowns, shoe covers, boots, respirators, face shields, safety glasses, or goggles. Individuals should be trained according to applicable regulatory and company/ institution requirements before working with potentially biohazardous materials. Follow all applicable local, state/provincial, and/or national regulations. The following references provide general guidelines when handling biological samples in laboratory environment.

 U.S. Department of Health and Human Services, Biosafety in Microbiological and Biomedical Laboratories (BMBL), 5th Edition, HHS Publication No. (CDC) 21-1112, Revised December 2009; found at:

https://www.cdc.gov/labs/pdf/CDC-BiosafetymicrobiologicalBiomedicalLaboratories-2009-P.pdf

 World Health Organization, Laboratory Biosafety Manual, 3rd Edition, WHO/CDS/CSR/LYO/2004.11; found at:

www.who.int/csr/resources/publications/biosafety/Biosafety7.pdf



Documentation and support

Related documentation

Document	Publication number
VeritiPro [™] Thermal Cycler PCR Starter Kit User Guide	MAN0019493
VeritiPro [™] Thermal Cycler Quick Reference (96-well)	MAN0019367
VeritiPro [™] Thermal Cycler Quick Reference (384-well)	MAN0024775

Customer and technical support

Visit thermofisher.com/support for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- Order and web support
- Product documentation
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

