

© Copyright 2007, 2010 Applied Biosystems. All rights reserved.

#### For Research Use Only. Not for use in diagnostic procedures.

Notice to Purchaser

The Applied Biosystems 7900HT Fast Real-Time PCR System is a real-time thermal cycler covered by US patents and corresponding claims in their non-US counterparts, owned by Applied Biosystems. No right is conveyed expressly, by implication or by estoppel under any other patent claim, such as claims to apparatus, reagents, kits, or methods such as 5' nuclease methods. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

#### TRADEMARKS:

ABI PRISM, Applied Biosystems, MicroAmp, and VIC are registered trademarks and AB (Design), Applera, FAM, JOE, NED, ROX, TAMRA, and TET are trademarks of Applied Biosystems or its subsidiaries in the U.S. and/or certain other countries.

GeneAmp and TaqMan are registered trademarks of Roche Molecular Systems, Inc.

SYBR is a registered trademark of Molecular Probes, Inc.

Zymark is a registered trademark of Zymark Corporation.

Windows is a registered trademark of Microsoft Corporation.

All other trademarks are the sole property of their respective owners.

Part Number 4365542 Rev. C 06/2010

# Contents

	Preface	V
	How to Use This Guide	v
Chapter 1	Performing Calibration and Verification Runs	1
	Recommended Run Schedule	
	Performing an Instrument Verification Run	32
Chapter 2	Maintaining the 7900HT Instrument Hardware	51
	Recommended Maintenance Schedule Removing and Installing a Sample Block Replacing the Plate Adapter Decontaminating the Sample Block	53 62
Chapter 3	Maintaining the Computer and Software	71
	Recommended Maintenance Schedule	
	Archiving and Backing Up SDS Software Files	
	Managing Local User Accounts in the SDS Software	78
	Updating the Operating System Software	
	Updating the Operating System Software	
Chapter 4	Maintaining the Automation Accessory	87
	Recommended Maintenance Schedule	
	Automation Accessory Components and Stack Positions	
	Adjusting the Plate Sensor Switch	
	Aligning the Plate Handler	97

	Aligning the Fixed-Position Bar Code Reader	
Chapter 5	Troubleshooting Chemistry and Assay Runs	119
	Low Precision or Irreproducibility	3 124
Chapter 6	Flags and Filtering	127
	Using Flags	128
	Flags By Analysis	
	Description of Flags	132
Appendix A	Adding Custom Dyes to the Pure Dye Set	141
	Before You Begin	141
	Creating a Dilution Series Plate	
	Creating a Plate Document for the Dilution Series Plate	
	Running and Analyzing the Dilution Series Plate	
	Creating a Pure Dye Plate with the Custom Dye(s)	
	Creating a Plate Document Template with the Custom Dye(s)	
Appendix B	Instrument Connections	153
Appendix C	Parts and Consumables	157
	Interchangeable Sample Block Modules and Accessories	157
	Consumables and Disposables	
	Instrument Maintenance and Verification	160
	Index	163

#### **Preface**

#### How to Use This Guide

### Purpose of This Guide

This guide provides recommended schedules and procedures for maintaining the Applied Biosystems 7900HT Fast Real-Time PCR System (7900HT Fast System). This guide also includes troubleshooting information for the maintenance procedures.

#### Audience

This guide is intended for:

- Personnel who maintain the 7900HT Fast System hardware
- System Administrators who maintain the 7900HT Fast System software
- Personnel running experiments

#### **Assumptions**

For hardware maintenance, this guide assumes that you have:

- Familiarity with the Microsoft Windows® XP operating system.
- Knowledge of general techniques for handling DNA samples and preparing them for PCR.

For software maintenance, this guide assumes that you have:

- Familiarity with the Microsoft Windows® XP operating system.
- A general understanding of hard drives and data storage, file transfers, and copying and pasting.
- Networking experience, if you want to integrate the 7900HT Fast System into your existing laboratory data flow

#### **Text Conventions**

This guide uses the following conventions:

- **Bold** indicates user action. For example:
  - Type 0, then press Enter for each of the remaining fields.
- *Italic* text indicates new or important words and is also used for emphasis. For example:
  - Before analyzing, always prepare fresh matrix.
- A right arrow bracket (>) separates successive commands you select from a dropdown or shortcut menu. For example:

Select File > Open.

## User Attention Words

The following user attention words appear in Applied Biosystems user documentation. Each word implies a particular level of observation or action as described below:

**Note** – Provides information that may be of interest or help but is not critical to the use of the product.

**IMPORTANT!** – Provides information that is necessary for proper instrument operation, accurate chemistry kit use, or safe use of a chemical.

**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.

#### Safety

Follow specific safety practices when using this instrument. For safety guidelines, refer to the "Safety and EMC Compliance" section in the *Applied Biosystems 7900HT Fast Real-Time PCR System Site Preparation Guide* (PN 4351923).

For any chemical manufactured or distributed by Applied Biosystems, you can obtain the MSDS from Applied Biosystems. This service is available free 24 hours a day. To obtain MSDSs:

- 1. Go to www.appliedbiosystems.com, click Support, then click MSDS Search.
- 2. In the Keyword Search field, enter the chemical name, product name, MSDS part number, or other information that appears in the MSDS of interest. Select the language of your choice, then click **Search**.
- 3. Find the document of interest, right-click the document title, then select any of the following:
  - **Open** To view the document
  - **Print Target** To print the document
  - Save Target As To download a PDF version of the document to a destination that you choose

For the MSDSs of chemicals not distributed by Applied Biosystems, contact the chemical manufacturer.

#### **How to Obtain More Information**

### Related Documentation

For more information about using the 7900HT Fast System, refer to:

- Sequence Detection Systems Software version 2.3 Online Help (SDS Online Help)
- Applied Biosystems 7900HT Fast Real-Time PCR System Site Preparation Guide (PN 4351923)
- Applied Biosystems 7900HT Fast Real-Time PCR System Allelic Discrimination Getting Started Guide (PN 4364015)
- Applied Biosystems 7900HT Fast Real-Time PCR System Absolute Quantitation Using Standard Curve Getting Started Guide (PN 4364014)
- Applied Biosystems 7900HT Fast Real-Time PCR System Plus-Minus Getting Started Guide (PN 4364017)
- Applied Biosystems 7900HT Fast Real-Time PCR System Relative Quantitation Using Comparative C<sub>T</sub> Getting Started Guide (PN 4364016)
- Real-Time PCR Systems Chemistry Guide (PN 4348358)

- SDS Enterprise Database for the Applied Biosystems 7900HT Fast Real-Time PCR System Administrators Guide (PN 4351669)
- TaqMan® Low Density Array Getting Started Guide (PN 4319399)

#### Send Us Your Comments

Applied Biosystems welcomes your comments and suggestions for improving its user documents. You can e-mail your comments to:

techpubs@appliedbiosystems.com

#### **How to Obtain Support**

To contact Applied Biosystems Technical Support from North America by telephone, call **1.800.762.4001**.

For the latest services and support information for all locations, go to <a href="http://www.appliedbiosystems.com">http://www.appliedbiosystems.com</a>, then click the link for Support.

At the Support page, you can:

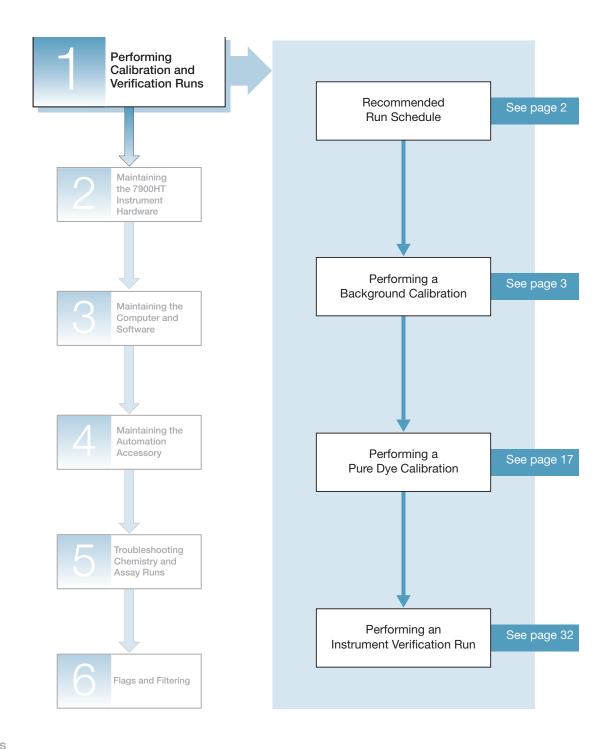
- Obtain worldwide telephone and fax numbers to contact Applied Biosystems Technical Support and Sales facilities
- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support
- Order Applied Biosystems user documents, MSDSs, certificates of analysis, and other related documents
- · Download PDF documents
- Obtain information about customer training
- Download software updates and patches

#### Preface

How to Obtain Support

# 1

# Performing Calibration and Verification Runs



#### Recommended Run Schedule

To ensure optimal performance of your 7900HT Fast System, perform the calibration and verification runs as indicated in the table below.

Schedule	Run Type
Semi-annually	Perform a background calibration (see page 3)
January July SuM T WTh F S SUM T WTh F S	<b>IMPORTANT!</b> You must also perform a background calibration before performing a pure dye calibration and after installing a new, uncalibrated sample block.
6 Months	Perform a pure dye calibration (see page 17)  IMPORTANT! You must also perform a pure dye calibration before performing an instrument verification run and after installing a new, uncalibrated sample block.
	Perform an instrument verification run (see page 32)

#### **Performing a Background Calibration**

#### Overview

#### When to Perform

Applied Biosystems recommends performing a background calibration:

- Every 6 months, or as often as needed depending on instrument use
- Before performing a pure dye calibration (see page 17)
- After installing a new, uncalibrated sample block (see page 53)

You can also perform background calibrations to detect and troubleshoot sample block contamination.

# About the Background Calibration

A background calibration measures the level of ambient fluorescence in the 7900HT instrument. During the run, the 7900HT instrument performs a continuous scan of a background plate (which contains PCR buffer or deionized water) for 2 minutes at 60 °C. Afterwards, the SDS software averages the spectrum recorded during the run and extracts the resulting spectral component to a calibration file. The SDS software uses the calibration file during subsequent runs to remove the background signal from the run data.

# About the Background Component

Fluorescence collected by the 7900HT Fast System includes a fluorescent signal inherent to the system, commonly referred to as background fluorescence. This background component is a composite signal found in all spectral data that consists of fluorescence from several sources, including: the background electronic signal, the sample block, the water inside the consumable, and the consumable itself. Because the background fluorescence can interfere with the precision of the data collected by the 7900HT System, the instrument is engineered to minimize the background signal; additionally, the SDS software algorithmically eliminates the background signal from each fluorescent sample to maximize the instrument's sensitivity.

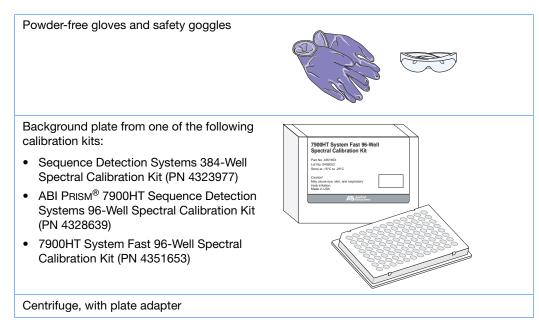
Notes\_\_\_\_\_

#### Before You Begin

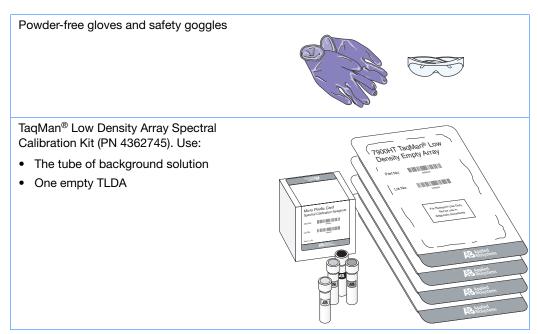
Time Required 30 minutes

Materials Required You need the materials listed below to perform this procedure.

#### For 384- or 96-Well Sample Blocks



#### For the TaqMan Low Density Array Block



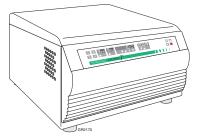
#### Vortex

Micropipettors and pipet tips:

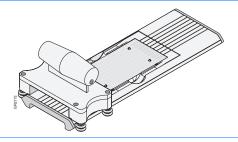
- Rainin F-100 micropipette (100-μL)
- Rainin Fine Point pipet tips (100-μL)



Centrifuge (For more information, see the TaqMan® Low Density Array Getting Started Guide.)



TaqMan<sup>®</sup> Low Density Array Sealer (For more information, see the *TaqMan*<sup>®</sup> Low Density Array Getting Started Guide.)



#### Workflow

#### To perform the background calibration:

- **1.** Obtain the tools and materials required (see pages 4 and 5).
- **2.** Prepare the background plate (see page 6) or background TLDA (see page 9).
- **3.** Create a plate document (see page 10).
- **4.** Run the background plate or background TLDA (see page 12).
- **5.** Analyze the background data (see page 14).

#### **Troubleshooting**

If you experience problems with the background calibration, see "Troubleshooting Background Calibrations" on page 15.

#### Preparing the Background Plate

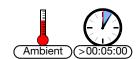
**IMPORTANT!** Wear powder-free gloves when you handle the background plate.



- **1.** Retrieve the calibration kit from the freezer and remove the packaged background plate from within it.
- **2.** Return the calibration kit to the freezer.



**3.** Allow the background plate to warm to room temperature (approximately 5 min).



**4.** Remove the background plate from its packaging.

**IMPORTANT!** Do not discard the packaging for the background plate. The background plate can be used repeatedly if it is stored in its original packaging sleeve.



**5.** Briefly centrifuge the background plate in a centrifuge with plate adapter.



	0			

**6.** Verify that the liquid in each well of the background plate is positioned at the bottom of the well. If not, centrifuge the plate again at a higher rpm and for a longer period of time.

**IMPORTANT!** Do not allow the bottom of the background plate to become dirty. Fluids and other contaminants that adhere to the bottom of the plate can contaminate the sample block and cause an abnormally high background signal.



Correct	Incorrect		
Sample is positioned at bottom of well.	The plate has not been centrifuged with enough force, or has not been centrifuged for enough time.		

**7.** Proceed to "Creating a Background Plate Document" on page 10.

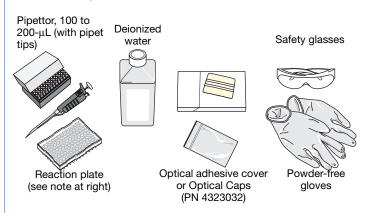
N	Л	$\overline{}$	+	$\overline{}$	0
ı,	И	U	U	U	J



#### Creating a Background Plate

If a background plate from a calibration kit is not available, you can create one to perform the background calibration by following the procedure below. Whenever possible, Applied Biosystems recommends the use of the background plate included with the calibration kit because it contains a PCR buffer that more accurately models the reagents used for PCR, and therefore produces better calibration data.

#### **Materials Required**



Note: Reaction plate options are:

- 384-Well Clear Optical Reaction Plate (standard 384-well reaction plate)
- MicroAmp® 96-Well Optical Reaction Plate (standard 96-well reaction plate)
- Optical 96-Well Fast Thermal Cycling Plate (Fast 96-well reaction plate)

#### To create the background plate:

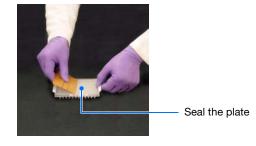
**IMPORTANT!** Wear powder-free gloves while creating the background plate.



- 1. Remove a reaction plate from its box and place it on a clean, dry
- Aliquot deionized water to each well of the reaction plate, as follows:
  - $-\,$  20  $\mu L$  per well for a standard 384-well or Fast 96-well reaction plate
  - 50 μL per well for a standard 96-well reaction plate

Pipet deionized water to each well

- 3. Seal the plate:
  - For standard 384-well or Fast 96-well reaction plates, use only an optical adhesive cover.
  - For standard 96-well reaction plates, use an optical adhesive cover or optical flat caps.



4. Continue with step 5 on page 6.

ΝI	_	ā.	_	_
IΜ	( )	ш	ш	5

#### Preparing the Background TLDA

- **1.** Retrieve the calibration kit from the freezer and remove:
  - The tube of background solution
  - · One empty TLDA
- **2.** Return the calibration kit to the freezer.



**3.** Allow the background solution to thaw at room temperature.



- **4.** When the solution has thawed, vortex the tube.
- 5. Load the background solution into the empty TLDA, loading 100  $\mu$ L of solution per fill reservoir.

**Note:** For detailed loading procedures see the *TaqMan® Low Density Array Getting Started Guide*.

**6.** Centrifuge and seal the background TLDA.

**Note:** For centrifuging and sealing procedures see the *TaqMan*<sup>®</sup> *Low Density Array Getting Started Guide*.

**7.** Proceed to "Creating a Background Plate Document" on page 10.



#### Creating a Background Plate Document

- Double-click on the desktop (or select Start > All Programs > Applied Biosystems > SDS
   2.3 > SDS
   2.3 > software.
- 2. If the login option is enabled, the Login dialog box appears. Enter your **User Name** and **Password**, then click **OK**.

**Note:** If the login option is not enabled, no Login dialog box appears. Skip to step 3 below.

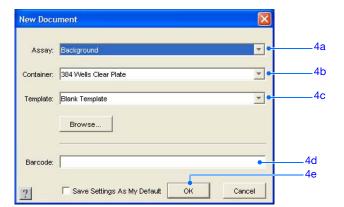


- 3. Click □ (or select File > New).
- **4.** Complete the New Document dialog box:
  - a. Assay Select Background.
  - **b.** Container Select the appropriate format.
  - **c.** Template Select the appropriate template, as described in the table below:

For Container	Select Template
384 Wells Clear Plate	Blank Template
96 Wells Clear Plate (for a Standard 96-Well Sample Block)	Blank Template
96 Wells Clear Plate (for a Fast 96-Well Sample Block)	Fast 96 Well Background Plate.sdt
384 Wells TaqMan Low Density Array	TaqMan Low Density Array Background Plate.sdt

- d. If the background plate or background TLDA is labeled with a barcode, click the **Barcode** field, then type in or scan the barcode number.
- e. Click or leave of the software creates and opens a plate document with the attributes for a background calibration.

**IMPORTANT!** Do not modify the background plate document. The method for a background calibration is coded into the SDS software and consists of a single

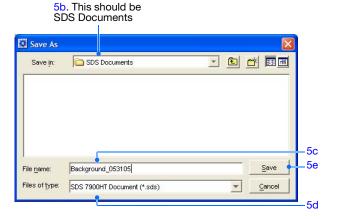


hold at 60 °C for 2 min. Because the plate contains only PCR buffer or deionized water, the plate document does not require sample or detector labels.

- **5.** Save the plate document:
  - a. Click (or select File ▶ Save) to open the Save As dialog box.
  - b. If the Save in field does not display SDS
     Documents, navigate to Applied
     Biosystems ▶ SDS Documents.
  - **c.** In the File name field, type in an appropriate file name, as described in the table below.

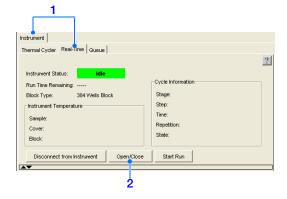
For Container	Туре
384 Wells Clear Plate	Background_ <date format="" in="" mmddyy=""></date>
	For example, the file name for a plate run on May 31, 2005, would be: <b>Background_053105</b> .
96 Wells Clear Plate (for a Standard 96-Well	Background_ <date format="" in="" mmddyy=""></date>
Sample Block)	For example, the file name for a plat run on May 31, 2005, would be: <b>Background_053105</b> .
96 Wells Clear Plate (for a Fast 96-Well Sample	FastBackground_Plate_ <date in<br="">MMDDYY format&gt;</date>
Block)	For example, the file name for a Fast plate run on May 31, 2005, would be: FastBackground_053105.
384 Wells TaqMan Low Density Array	Background <date format="" in="" mmddyy="">_TLDA</date>
	For example, the file name for a background TLDA run on May 31, 2005, would be:  Background_053105_TLDA.

- d. Select **SDS 7900HT Document (\*.sds)** from the Files of type drop-down list.
- e. Click \_\_\_\_\_\_. The software saves the plate document.
- **6.** Continue with "Running the Background Plate or Background TLDA" on page 12.



#### Running the Background Plate or Background TLDA

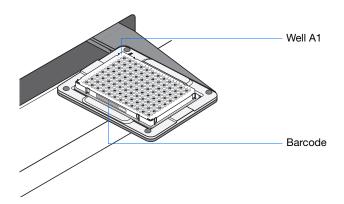
- 1. In the background plate document, select the **Instrument** > **Real-Time** tabs.
- **2.** Click Open/Close The instrument tray rotates to the OUT position.



**3.** Place the background plate or background TLDA into the instrument tray as shown.

**IMPORTANT!** The background plate or TLDA must be oriented so that the A1 position matches the A1 label on the instrument tray.

Standard 384-well reaction plate



**4.** Click Start Run. The instrument tray rotates to the IN position and performs the background calibration.

**Note:** Before starting the run, the instrument may pause (up to 15 min) to bring the heated cover to the appropriate temperature.

5. When the background calibration is complete and the Run Complete dialog box appears, click or lose the dialog box.

**6.** Click open/Close , then remove the background plate or background TLDA from the instrument tray.

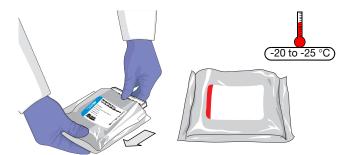
WARNING PHYSICAL INJURY

**HAZARD**. During instrument operation, the sample block can be heated as high as 100 °C. Before removing the background plate or background TLDA, be sure to wait until the sample block reaches room temperature.

**7.** Place the background plate or background TLDA back inside its packaging, then return it to the calibration kit in the freezer.

**IMPORTANT!** Do not discard the background plate or background TLDA. If you store it in its original packaging, you can use the background plate or background TLDA for up to 6 months after you open it.

**8.** Continue with "Analyzing the Background Data" on page 14.

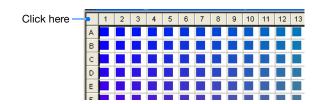


#### Analyzing the Background Data

When you analyze the background data, the SDS software extracts the calibration values from the background plate document, then provides pass/fail information. After extraction, the SDS software stores the data as part of the calibration file located in the Calibration subdirectory of the SDS directory.

#### To analyze the background data:

1. In the Plate Grid of the background plate document, click the box above the A label to select all wells.



#### 2. Select Analysis > Extract Background.

The software attempts to extract the background signal and displays the success of the extraction in a dialog box.

**3.** If the software displays:

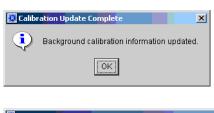
Calibration Update Complete – The analysis is successful. The raw spectra read from the background plate conform to acceptable limits.

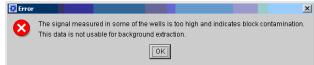
Click \_\_\_\_, then go to step 4.

*Error* – The run is unsuccessful. The software stopped the extraction because one or more raw spectra exceed 4000 FSU.

Click \_\_\_\_\_, then troubleshoot the failed run as explained in "Troubleshooting Background Calibrations" on page 15.

- **4.** Click (or select **File ▶ Save**) to save the background plate document.
- **5.** Select **File** ▶ **Close** to close the background plate document.





#### **Troubleshooting Background Calibrations**

### Troubleshooting Table

Table 1-1 Troubleshooting background calibrations

Observation	Possible Cause	Recommended Action	
Software will not extract background data	During setup, the wrong plate type was assigned to the plate document	Run a new background plate document with the proper plate type setting.	
	Background is too high (>4000 FSU <sup>‡</sup> )	See "Isolating Sample Block Contamination" below.	
Background is too high (>4000 FSU)	Sample block contamination	Construct and run a new background plate.	
	Background plate contamination	2. See "Isolating Sample Block Contamination" below.	

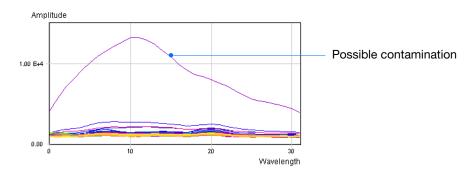
<sup>‡</sup>Fluorescent standard units – The measure of amplitude displayed along the Y-axis of the Background Plot.

#### Isolating Sample Block Contamination

Signals exceeding 4000 FSU are considered outside the limit of normal background fluorescence and indicates that the either the background plate or the sample block may be contaminated.

- 1. Open the plate document for the background calibration.
- 2. In the toolbar, click [ (Hide/Show System Raw Data Pane).
- Select all wells in the plate document.
   The SDS software highlights the selected wells and displays the raw spectral data.
- 4. Inspect the raw background data for an aberrant spectral peak or peaks.

  Wells producing raw spectra that exceed 4000 FSU are considered irregular and could be contaminated. The following figure illustrates the raw data produced by a run on a sample block module containing a contaminated well.



**5.** Drag the vertical bar in the temperature plot to inspect all the data points. The location(s) of the contaminated well(s) is displayed in tooltip when you move the mouse over the curves in the graph.

# **Chapter 1** Performing Calibration and Verification Runs Performing a Background Calibration



- 6. Decontaminate the sample block as explained on page 65.
- 7. Run a background plate to confirm that the contaminants have been removed.

  If the contamination is present after running the background plate for a second time, the background plate is likely to be the source of contamination.

Notes\_\_\_\_\_

#### Performing a Pure Dye Calibration

#### Overview

#### When to Perform

Applied Biosystems recommends performing a pure dye calibration:

• Every 6 months, depending on instrument use

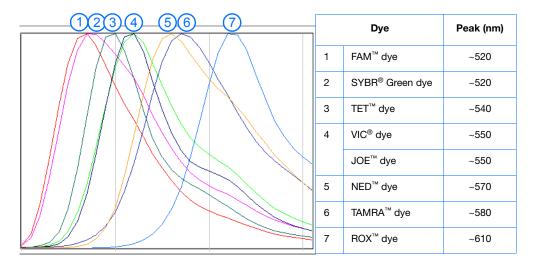
**IMPORTANT!** The age and use of instrument components can affect pure spectra readings. Update the pure spectra data files once or twice annually, depending on the frequency of instrument use.

- Before performing an instrument verification run (see page 32)
- After installing a new, uncalibrated sample block (see page 53)

#### Purpose of the Pure Dye Calibrations

Before operating the instrument, you must generate spectral data by performing a pure dye calibration. During the pure dye calibration, the SDS software collects spectral data from a set of pure dye standards during a 2-minute hold at 60 °C. The SDS software stores the spectral data in the pure spectra run file, a calibration file located in the SDS directory. The SDS software uses the spectral data from the set of pure dye standards to distinguish the individual contribution of each dye in the collective fluorescence gathered by the instrument during a run.

After each run, the SDS software receives run data in the form of a raw spectra signal for each reading. To make sense of the raw data, the software must determine the contribution of each fluorescent dye used in the sample by comparing the raw spectra to the spectral data contained in the pure spectra run file.



**Note:** When a plate document is saved after analysis, the SDS software stores the pure spectra information with the collected fluorescent data for that experiment.

# Chapter 1 Performing Calibration and Verification Runs Performing a Pure Dye Calibration

#### Components of the Pure Dye Standards

The 7900HT Fast System is calibrated with several dyes, which are included in the calibration kits.

- The Sequence Detection Systems 384-Well Spectral Calibration Kit is provided for the standard 384-well sample block. This kit includes one 384-well pure dye plate, preloaded with the following dyes: FAM<sup>™</sup> dye, JOE<sup>™</sup> dye, NED<sup>™</sup> dye, ROX<sup>™</sup> dye, SYBR<sup>®</sup> Green dye, TAMRA<sup>™</sup> dye, TET<sup>™</sup> dye, and VIC<sup>®</sup> dye.
- The ABI PRISM<sup>®</sup> 7900HT Sequence Detection Systems 96-Well Spectral Calibration Kit is provided for the standard 96-well sample block. This kit includes two 96-well pure dye plates. Plate 1 is preloaded with the following dyes: FAM<sup>™</sup> dye, JOE<sup>™</sup> dye, NED<sup>™</sup> dye, and ROX<sup>™</sup> dye. Plate 2 is preloaded with the following dyes: SYBR<sup>®</sup> Green dye, TAMRA<sup>™</sup> dye, TET<sup>™</sup> dye, and VIC<sup>®</sup> dye.
- The 7900HT System Fast 96-Well Spectral Calibration Kit is provided for the Fast 96-well sample block. This kit includes two 96-well pure dye plates. Plate 1 is preloaded with the following dyes: FAM<sup>™</sup> dye, JOE<sup>™</sup> dye, NED<sup>™</sup> dye, and ROX<sup>™</sup> dye. Plate 2 is preloaded with the following dyes: SYBR<sup>®</sup> Green dye, TAMRA<sup>™</sup> dye, TET<sup>™</sup> dye, and VIC<sup>®</sup> dye.
- The TaqMan<sup>®</sup> Low Density Array Spectral Calibration Kit is provided for the TaqMan<sup>®</sup> Low Density Array block. The kit includes three tubes of dye: FAM<sup>™</sup> dye, ROX<sup>™</sup> dye, and VIC<sup>®</sup> dye and three empty TLDAs. To perform a pure dye calibration, you load each dye into an empty TLDA.

#### Custom Pure Dyes

The 7900HT Fast System supports the detection of custom pure dyes (dyes other than those provided by Applied Biosystems). To add custom pure dyes to the pure dye set for your instrument, see Appendix A on page 141.

Notes		

#### **Before You Begin**

#### Time Required 30 minutes

#### Materials Required

You need the materials listed below to perform this procedure.

#### For 384- or 96-Well Sample Blocks

Pure dye plate(s) from one of the following calibration kits:

Sequence Detection Systems 384-Well Spectral Calibration Kit (PN 4323977)

ABI PRISM® 7900HT Sequence Detection Systems 96-Well Spectral Calibration Kit (PN 4328639)

7900HT System Fast 96-Well Spectral Calibration Kit (PN 4351653)

Note: Both the standard 96-well and Fast 96-well calibration kits contain two pure dye plates.

Centrifuge, with plate adapter

#### For the TaqMan Low Density Array Block

TaqMan® Low Density Array Spectral
Calibration Kit (PN 4362745). Use:

• The three dye tubes: FAM™ dye,
ROX™ dye, and VIC® dye

• Three empty TLDAs



#### Vortex

Micropipettors and pipet tips:

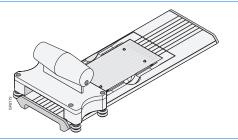
- Rainin F-100 micropipette (100-μL)
- Rainin Fine Point pipet tips (100-μL)



Centrifuge (For more information, see the *TaqMan*<sup>®</sup> *Low Density Array Getting Started Guide.*)



TaqMan® Low Density Array Sealer (For more information, see the *TaqMan® Low Density Array Getting Started Guide.*)



#### Workflow To perform the pure dye calibration:

**1.** Perform a background calibration first (see page 3)

**IMPORTANT!** You must perform a background calibration before every pure dye calibration.

- **2.** Obtain the tools and materials required (see pages 19 and 20).
- **3.** Prepare the pure dye plate(s) (see page 21) or pure dye TLDAs (see page 23).
- **4.** Create a plate document (see page 24).
- **5.** Run the pure dye plate(s) or pure dye TLDAs (see page 27).
- **6.** Analyze the pure dye data (see page 29).

#### **Troubleshooting**

If you experience problems with the pure dye calibration, see "Troubleshooting Pure Dye Calibrations" on page 31.

B. II				
I/I	$\sim$	·T	$\sim$	c

#### Preparing the Pure Dye Plate(s)

**IMPORTANT!** You must perform a background calibration before every pure dye calibration.

**IMPORTANT!** Wear powder-free gloves when you handle the pure dye plates.

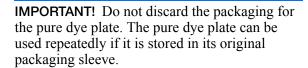


**1.** Retrieve the calibration kit from the freezer, then remove the packaged pure dye plate(s).

**Note:** The standard 384-well calibration kit contains one pure dye plate. Both the standard 96-well and Fast 96-well calibration kits contain two pure dye plates.

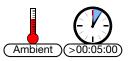
- 2. Return the calibration kit to the freezer.
- **3.** Allow the pure dye plate(s) to warm to room temperature (approximately 5 min).
- **4.** Remove a pure dye plate from its packaging.

**IMPORTANT!** Do not remove a pure dye plate from its packaging until you are ready to run it. The fluorescent dye contained in the wells of each pure dye plate is photosensitive. Prolonged exposure to light can diminish the fluorescent signal strength of the plate.



**5.** Briefly centrifuge the pure dye plate in a centrifuge with plate adapter.











**6.** Verify that the pure dye standard in each well of the pure dye plate is positioned at the bottom of the well. If not, centrifuge the plate again at a higher rpm and for a longer period of time.



Correct	Incorrect		
Sample is positioned at bottom of well.	The plate has not been centrifuged with enough force, or has not been centrifuged for enough time.		

**7.** Proceed to "Creating a Pure Dye Plate Document" on page 24.

#### Preparing the Pure Dye TLDAs

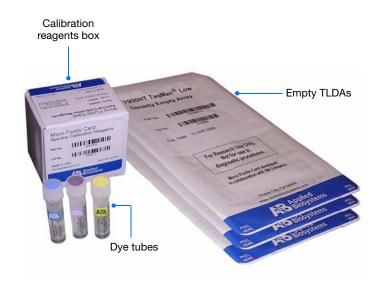
**IMPORTANT!** You must perform a background calibration before every pure dye calibration.

**IMPORTANT!** Wear powder-free gloves when performing this procedure.

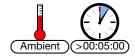


WARNING Spectral Calibration Bulk FAM<sup>TM</sup>, ROX<sup>TM</sup>, and VIC<sup>®</sup> Dyes are combustible liquids and vapors. Exposure causes eye, skin, and respiratory tract irritation. Read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

- **1.** Retrieve the calibration kit from the freezer and remove:
  - The three dye tubes: FAM dye, ROX dye, and VIC dye
  - Three empty TLDAs
- **2.** Return the calibration kit to the freezer.



**3.** Allow the dyes to thaw at room temperature.



- **4.** When the dyes have thawed, vortex the tubes.
- **5.** Load the FAM dye into one of the empty TLDAs, loading 100 μL of dye per fill reservoir.

**Note:** For detailed loading procedures see the *TaqMan® Low Density Array Getting Started Guide*.





- **6.** Repeat step 5 for the ROX and VIC dyes, using the two remaining empty TLDAs.
- **7.** Centrifuge and seal the pure dye TLDAs.

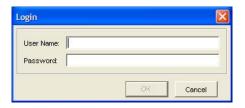
**Note:** For centrifuging and sealing procedures see the *TaqMan*<sup>®</sup> *Low Density Array Getting Started Guide*.

**8.** Proceed to "Creating a Pure Dye Plate Document" on page 24.

#### Creating a Pure Dye Plate Document

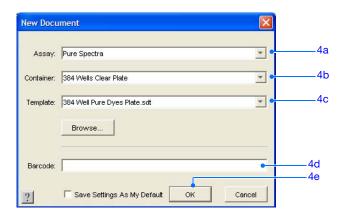
- Double-click on the desktop (or select Start > All Programs > Applied Biosystems > SDS
   2.3 > SDS
   2.3) to start the SDS software.
- 2. If the login option is enabled, the Login dialog box appears. Enter your User Name and Password, then click OK.

**Note:** If the login option is not enabled, no Login dialog box appears. Skip to step 3 below.



- 3. Click □ (or select File > New).
- **4.** Complete the New Document dialog box:
  - a. Assay Select Pure Spectra.
  - **b.** Container Select the appropriate format.
  - **c.** Template Select the appropriate template, as described in the table below.

For Container	Select Template		
384 Wells Clear Plate	384 Well Pure Dyes Plate.sdt		
96 Wells Clear Plate (for a Standard 96-Well Sample Block)	96 Well Pure Dyes Plate 1.sdt to run Plate 1 (containing FAM, JOE, NED, and ROX dyes)     96 Well Pure Dyes Plate 2.sdt to run Plate 2 (containing SYBR Green, TAMRA, TET, and VIC dyes)		



For Container	Select Template
96 Wells Clear Plate (for a Fast 96-Well Sample Block)	Fast 96 Well Pure Dyes Plate 1.sdt to run Plate 1 (containing FAM, JOE, NED, and ROX dyes)     Fast 96 Well Pure Dyes Plate 2.sdt to run Plate 2 (containing SYBR Green, TAMRA, TET, and VIC dyes)
384 Wells TaqMan Low Density Array	TaqMan Low Density Array     FAM Pure Dye.sdt to run the pure     dye TLDA containing FAM dye
	TaqMan Low Density Array     ROX Pure Dye.sdt to run the pure     dye TLDA containing ROX dye
	TaqMan Low Density Array     VIC Pure Dye.sdt to run the pure     dye TLDA containing VIC dye

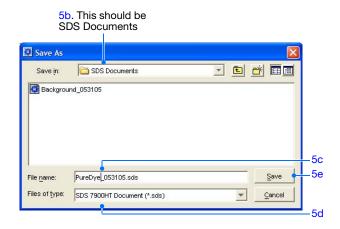
- d. If the pure dye plate or pure dye TLDA is labeled with a barcode, click the **Barcode** field, then type in or scan the barcode number.

**IMPORTANT!** Do not modify the pure dye plate document.

#### **5.** Save the plate document:

- a. Click ☐ (or select File ▶ Save) to open the Save As dialog box.
- b. If the Save in field does not display SDS
   Documents, navigate to Applied
   Biosystems > SDS Documents.
- **c.** In the File name field, type in an appropriate file name, as described in the table below.

For Container	Туре
384 Wells Clear Plate	PureDye_ <date format="" in="" mmddyy=""></date>
	For example, the file name for a plate run on May 31, 2005, would be: PureDye_053105.
96 Wells Clear Plate (for a Standard 96-Well Sample Block)	PureDye_Plate <plate #="">_<date format="" in="" mmddyy=""></date></plate>
	For example, the file name for a Plate 1 run on May 31, 2005, would be: PureDye_Plate1_053105.





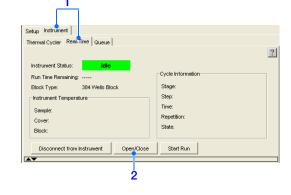
For Container	Туре	
96 Wells Clear Plate (for a Fast 96-Well Sample Block)	FastPureDye_Plate <plate #="">_<date format="" in="" mmddyy=""></date></plate>	
	For example, the file name for a Fast Plate 1 run on May 31, 2005, would be: FastPureDye_Plate1_053105.	
384 Wells TaqMan Low Density Array	<pre><dye><date format="" in="" mmddyy="">_TLDA</date></dye></pre>	
	For example, the file name for a FAM dye TLDA run on May 31, 2005, would be: <b>FAM_053105_TLDA</b> .	

- d. Select **SDS 7900HT Document (\*.sds)** from the Files of type drop-down list.
- **e.** Click \_\_\_\_\_\_. The software saves the plate document.
- **6.** Continue with "Running a Pure Dye Plate or Pure Dye TLDA" on page 27.

Notes			

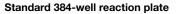
#### Running a Pure Dye Plate or Pure Dye TLDA

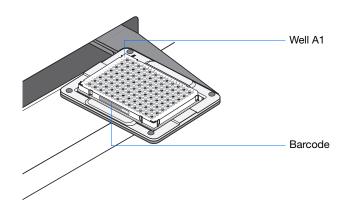
- 1. In the pure dye plate document, select the **Instrument** > **Real-Time** tabs.
- **2.** Click Open/Close The instrument tray rotates to the OUT position.



**3.** Place the pure dye plate or pure dye TLDA into the instrument tray as shown.

**IMPORTANT!** The pure dye plate or TLDA must be oriented so that the A1 position matches the A1 label on the instrument tray.





**4.** Click \_\_start Run\_. The instrument tray rotates to the IN position and performs the pure dye calibration.

**Note:** Before starting the run, the instrument may pause (up to 15 min) to bring the heated cover to the appropriate temperature.

5. When the pure dye calibration is complete and the Run Complete dialog box appears, click or to close the dialog box.





**6.** Click open/Close , then remove the pure dye plate or pure dye TLDA from the instrument tray.

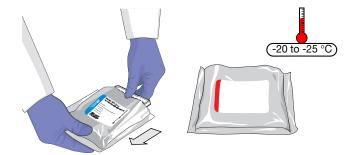


HAZARD. During instrument operation, the sample block can be heated as high as 100 °C. Before removing the background plate or background TLDA, be sure to wait until the sample block reaches room temperature.

**7.** Place the pure dye plate or each pure dye TLDA inside its packaging, then return it to the calibration kit in the freezer.

**IMPORTANT!** Do not discard the pure dye plate or pure dye TLDAs. If you store it in its original packaging, you can use the pure dye plate or each pure dye TLDA for up to 6 months after you open it.

**8.** Continue with "Analyzing the Pure Dye Data" on page 29.



# Analyzing the Pure Dye Data

When you analyze the pure dye data, the SDS software extracts the calibration values from the pure dye plate document, then displays the spectral data in the Pure Dye Wizard.

The purpose of viewing the data in the Pure Dye Wizard is to eliminate irregular pure dye peaks from the data set. The wizard presents the spectral data in sets of three wells, each containing the same pure dye. Because the wells displayed by the wizard contain the pure dye at an identical concentration, the signal peaks for the set should be identical. Occasionally, pipetting inaccuracies or contamination can cause a well signal to shift slightly. While viewing the data, you must eliminate any outlying peaks.

### To analyze the pure dye data:

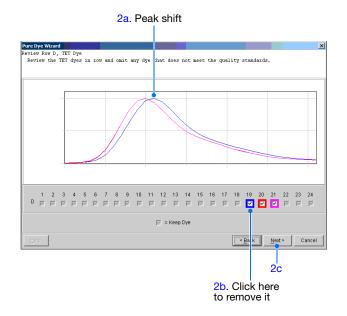
- 1. In the pure dye plate document, select Analysis ▶ Extract Pure Dye Wizard.
- **2.** Follow the instructions as explained in the Pure Dye Wizard to extract the pure dye spectra. When presented with each screen:
  - **a.** Inspect the spectra for shifts in peak location.
  - b. If the data set contains an outlying peak, eliminate it by clicking the check box of the associated well.

**Note:** Dye spectra are generally acceptable if they peak at the same location as their group but diverge slightly at other wavelengths.

- **c.** Click Next > 0 when finished.
- d. Repeat steps a through c for all remaining wells until prompted with a message reporting the extraction of the pure dyes.

The software extracts the pure spectra and stores the data as a component of the calibration file.

3. Click (or select File ▶ Save) to save the pure dye plate document.





- **4.** Select **File ▶ Close** to close the pure dye plate document.
- **5.** If you are performing the pure dye calibration for a:
  - Standard 384-well sample block The pure dye calibration is complete.
  - Standard 96-well or Fast 96-well sample block Run the second pure dye plate by repeating the following procedures:
    - "Creating a Pure Dye Plate Document" on page 24
    - "Running a Pure Dye Plate or Pure Dye TLDA" on page 27
    - "Analyzing the Pure Dye Data" on page 29
  - TaqMan Low Density Array Run the second and third pure dye TLDAs by repeating the following procedures for each TLDA:
    - "Creating a Pure Dye Plate Document" on page 24
    - "Running a Pure Dye Plate or Pure Dye TLDA" on page 27
    - "Analyzing the Pure Dye Data" on page 29

**IMPORTANT!** You must calibrate the instrument for all pure dye plates or pure dye TLDAs provided in your calibration kit.

Notes		

# **Troubleshooting Pure Dye Calibrations**

# Troubleshooting Table

Table 1-2 Troubleshooting pure dye calibrations

Observation	Possible Cause	Recommended Action				
Software will not extract pure dye data	During plate setup, the wrong plate type was assigned to the plate document	Create and run a new pure dye plate document with the proper plate type setting				
	A background plate was not run before the pure dye plate or TLDA	Run a background plate, then run the pure dye plate or TLDA again				
Raw data from pure dye calibration appears strange (see below)	Pure dye plate or TLDA was loaded backwards	Verify the pure dye     wavelengths are as     expected.				
		Rerun the pure dye plate or TLDA.				
Raw Data Plot   Calibration Data	<u>I</u>	<u>I</u>				
3.00 E44 2.50 E44 1.00 E44 5.00 E43 0.00 0 5 10 15 20 25 30						
2 59-1 Time						
Signals plateau (saturation) Signal is too low (<10,000 FSU)	Intensity is set too high/low	Call Applied Biosystems Technical Support.				
More than two outliers per dye in a single row	<ul><li>Evaporation</li><li>Contamination</li></ul>	Rerun the pure dye plate or TLDA. If the problem persists, discard the pure dye plate or TLDA and run a new one.				

# Performing an Instrument Verification Run

# Overview

### When to Perform

Applied Biosystems recommends performing an instrument verification run:

- Every 6 months
- As needed to verify the function of the 7900HT Fast System

# Purpose of the Verification Run

An instrument verification run is used to verify the performance of the 7900HT Fast System. During the run, TaqMan® RNase P reagents detect and quantify genomic copies of the human RNase P gene. (The RNase P gene is a single-copy gene encoding the RNA moiety of the RNase P enzyme.)

For the 7900HT Fast System instrument verification runs, the TaqMan RNase P reagents are available in the following formats:

- TaqMan® RNase P Instrument Verification Plates (RNase P plates) Verifies instrument performance for 7900HT instruments that use a Standard 384-Well Block, Standard 96-Well Block, or Fast 96-Well Block.
- TaqMan® Low Density Array RNase P Kit (TLDA RNase P kit) Verifies
  instrument performance for 7900HT instruments that use a TaqMan® Low Density
  Array Block.

# About RNase P Plates

RNase P plates are pre-loaded with the reagents required to detect and quantify genomic copies of the human RNase P gene. Each well contains pre-loaded reaction mix (TaqMan<sup>®</sup> Universal PCR Master Mix, RNase P primers, and FAM<sup>™</sup> dye-labeled probe) and a known concentration of human genomic DNA template.

Notes		

The table below illustrates the arrangement of standards and samples on each type of RNase P plate that is available for the 7900HT Fast System. The RNase P plates contain five replicate groups of standards (1250, 2500, 5000, 10,000, and 20,000 copies), two unknown populations (5000 and 10,000 copies), and template control (NTC) wells.

RNase P Plate	Sample Configuration
TaqMan <sup>®</sup> RNase P 384-Well Instrument Verification Plate	Unknown 1  5000  1
TaqMan <sup>®</sup> RNase P Instrument Verification Plate	1 2 3 4 5 6 7 8 9 10 11 12 A O O O O O O O O O O O O O O O O O O O
TaqMan <sup>®</sup> RNase P Fast 96- Well Instrument Verification Plate	B   5000   5000   STD 2500   E   STD 5000   STD 10000   STD 20000   STD 2000

# **Chapter 1** Performing Calibration and Verification Runs Performing an Instrument Verification Run



# About the TLDA RNase P Kit

The TLDA RNase P Kit includes one empty TLDA and eight tubes of solution. Each tube of solution contains reaction mix (TaqMan<sup>®</sup> Universal PCR Master Mix, RNase P primers, and FAM<sup>™</sup>-MGB dye-labeled probe) and a known concentration of human genomic DNA template.

To perform an instrument verification run, you load each solution into the empty TLDA. The figure below illustrates the arrangement of standards and samples. There are five replicate groups of standards (200, 400, 800, 1600, and 3200 copies), two unknown populations (800 and 1600 copies), and no template control (NTC).

	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	PORT
Α										N.	тС														1
В																									
С										Ш	nkr	101	vn	Δ (	ጸበ	U)									2
D	$ldsymbol{ld}}}}}}$									<u> </u>			•	,,,	00	٠,								Ш	
Е										H	nkr	101	vn	R	116	00	١								3
F										٠.			• • • •		,.,	-	_								
G										St	an	da	rd	200	<b>b</b> —										4
Н	$ldsymbol{ldsymbol{eta}}$												_		_										
ı										St	an	da	rd	400	<b>b</b> —										- 5
J	$oxed{oxed}$												_		_										
K										St	an	da	rd	800	<b>b</b> —										6
L	_									_			_	-	_										
M	_									St	an	da	rd	160	00										7
N	ㄴ												_		_										
0										St	an	da	rd	320	00										- 8
Р										-					Ţ <b>"</b>										

Notes\_\_\_\_\_



# **Before You Begin**

#### Time Required 2 hours

# Materials Required

You need the materials listed below to perform this procedure.

### For 384- or 96-Well Sample Blocks

Powder-free gloves and safety goggles Appropriate RNase P plate for your sample block: TaqMan® • TaqMan® RNase P 384-Well Instrument Verification Plate (PN 4323306) AB Applied Biosyste TaqMan® RNase P Instrument Verification Plate (PN 4310982, standard 96-well reaction plate) TagMan® RNase P Fast 96-Well Instrument Verification Plate (PN 4351979) Centrifuge, with plate adapter

# For the TaqMan Low Density Array Block

Powder-free gloves and safety goggles TaqMan® Low Density Array RNase P Kit (PN 4351468) 7900HT TaqMan® Low



### Micropipettors and pipet tips:

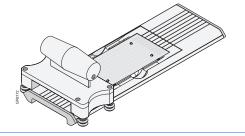
- Rainin F-100 micropipette (100-μL)
- Rainin Fine Point pipet tips (100-μL)



Centrifuge (For more information, see the *TaqMan® Low Density Array Getting Started Guide.*)



TaqMan<sup>®</sup> Low Density Array Sealer (For more information, see the *TaqMan*<sup>®</sup> Low Density Array Getting Started Guide.)



# Workflow To perform an instrument verification run:

- **1.** Perform a pure dye calibration (see page 17) before performing the instrument verification run.
- **2.** Obtain the tools and materials required (see pages 35 and 36).
- **3.** Prepare the RNase P plate (see page 37) or RNase P TLDA (see page 39).
- **4.** Create a plate document (see page 41).
- **5.** Run the RNase P plate or RNase P TLDA (see page 43).
- **6.** Analyze the run data (see page 44).
- 7. Verify instrument performance (see page 47).

# Preparing the RNase P Plate

**IMPORTANT!** Wear powder-free gloves when you handle the RNase plate.

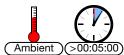


**1.** Retrieve the RNase P package from the freezer, then remove the packaged RNase P plate.

**Note:** For the standard 96-well RNase P plate, also use the compression pad provided in the package. Compression pads are not provided and should not be used for the Fast 96-well or standard 384-well RNase P plates.



**2.** Allow the RNase P plate to warm to room temperature (approximately 5 min).



**3.** Remove the RNase P plate from its packaging.



**4.** Briefly centrifuge the RNase P plate in a centrifuge with plate adapter.





**5.** Verify that the liquid in each well of the RNase P plate is positioned at the bottom of the well. If not, centrifuge the plate again at a higher rpm and for a longer period of time.



Correct	Incorrect
Sample is positioned at bottom of well.	The plate has not been centrifuged with enough force, or has not been centrifuged for enough time.

**6.** Proceed to "Creating a Plate Document for the Verification Run" on page 41.

**IMPORTANT!** To prevent degradation of the reaction components, start the run as soon as possible.

# Preparing the RNase P TLDA

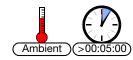
**IMPORTANT!** Wear powder-free gloves when performing this procedure.



- **1.** Remove the TLDA RNase P kit from the freezer, then unpack the contents:
  - Eight tubes of solution
  - · One empty TLDA
- **2.** Remove the empty TLDA from its packaging.



**3.** Allow the eight tubes of solution to thaw at room temperature (approximately 5 min).



- **4.** Mix the contents of each tube by tapping gently.
- **5.** Using the empty TLDA, load 100  $\mu$ L of the first solution into the designated reservoir.

**Note:** The insert in the TLDA RNase P kit describes the designated reservoirs.

**Note:** For detailed loading procedures see the *TaqMan® Low Density Array Getting Started Guide*.

**6.** Repeat step 5 for the remaining solutions.

# Chapter 1 Performing Calibration and Verification Runs Performing an Instrument Verification Run

**7.** Centrifuge and seal the three RNase P TLDAs.

**Note:** For centrifuging and sealing procedures see the *TaqMan*<sup>®</sup> *Low Density Array Getting Started Guide*.

**8.** Proceed to "Creating a Plate Document for the Verification Run" on page 41.

**IMPORTANT!** To prevent degradation of the reaction components, start the run as soon as possible.

Notes\_\_\_\_

# Creating a Plate Document for the Verification Run

- Double-click on the desktop (or select Start > All Programs > Applied Biosystems > SDS
   2.3 > SDS
   2.3) to start the SDS software.
- 2. If the login option is enabled, the Login dialog box appears. Enter your User Name and Password, then click OK.

**Note:** If the login option is not enabled, no Login dialog box appears. Skip to step 3 below.

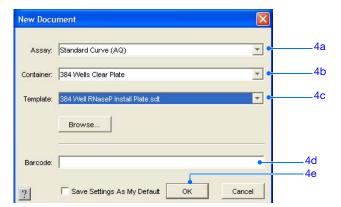


- 3. Click □ (or select File > New).
- **4.** Complete the New Document dialog box:
  - a. Assay Select Standard Curve (AQ).
  - **b.** Container Select the appropriate format.
  - **c.** Template Select the appropriate template, as described in the table below.

For Container	Select Template
384 Wells Clear Plate	384 Well RNaseP Install Plate.sdt
96 Wells Clear Plate (for a Standard 96-Well Sample Block)	96 Well RNaseP Install Plate.sdt
96 Wells Clear Plate (for a Fast 96-Well Sample Block)	Fast 96 Well RNaseP Install.sdt
384 Wells TaqMan Low Density Array	TaqMan Low Density Array RNaseP.sdt

- d. If the RNase P plate or RNase P TLDA is labeled with a barcode, click the **Barcode** field, then type in or scan the barcode number.
- e. Click \_\_\_ok\_\_\_. The software creates and opens a plate document.

**IMPORTANT!** Do not modify the plate document. The plate document template is programmed with the appropriate detector and method information for the instrument verification run.



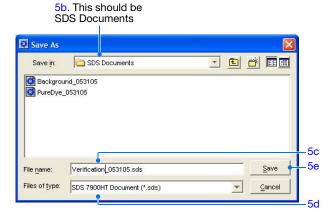
# **Chapter 1** Performing Calibration and Verification Runs Performing an Instrument Verification Run



- **5.** Save the plate document:
  - a. Click (or select File ▶ Save) to open the Save As dialog box.
  - b. If the Save in field does not display SDS
     Documents, navigate to Applied
     Biosystems ▶ SDS Documents.
  - **c.** In the File name field, type in an appropriate file name, as described in the table below.

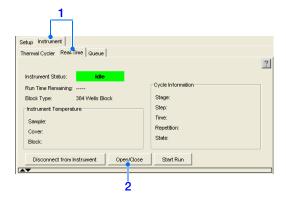
For Container	Type
384 Wells Clear Plate	Verification_ <date format="" in="" mmddyy=""></date>
	For example, the file name for a plate run on May 31, 2005, would be: <b>Verification_053105</b> .
96 Wells Clear Plate (for a Standard 96-Well	Verification_ <date format="" in="" mmddyy=""></date>
Sample Block)	For example, the file name for a plate run on May 31, 2005, would be: <b>Verification_053105</b> .
96 Wells Clear Plate (for a Fast 96-Well Sample	FastVerification_ <date format="" in="" mmddyy=""></date>
Block)	For example, the file name for a Fast plate run on May 31, 2005, would be: FastVerifiation_053105.
384 Wells TaqMan Low Density Array	Verification <date format="" in="" mmddyy="">_TLDA</date>
	For example, the file name for an RNase P TLDA run on May 31, 2005, would be: <b>Verification_053105_TLDA</b> .

- d. Select **SDS 7900HT Document (\*.sds)** from the Files of type drop-down list.
- e. Click \_\_save \_\_. The software saves the plate document.
- **6.** Continue with "Running the RNase P Plate or RNase P TLDA" on page 43.



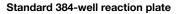
# Running the RNase P Plate or RNase P TLDA

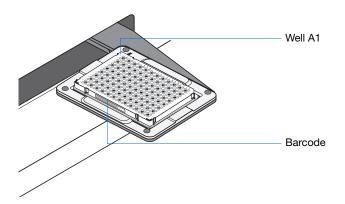
- 1. In the plate document, select the **Instrument** ▶ **Real-Time** tabs.
- **2.** Click Open/Close The instrument tray rotates to the OUT position.



**3.** Place the RNase P plate or RNase P TLDA into the instrument tray as shown.

**IMPORTANT!** The RNase P plate or TLDA must be oriented so that the A1 position matches the A1 label on the instrument tray.





**4.** Click Start Run. The instrument tray rotates to the IN position and performs the instrument verification run.

**Note:** Before starting the run, the instrument may pause (up to 15 min) to bring the heated cover to the appropriate temperature.

**5.** When the instrument verification run is complete and the Run Complete dialog box appears, click to close the dialog box.



**6.** Click open/Close , then remove the RNase P plate or RNase P TLDA from the instrument tray.

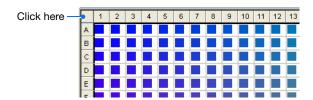


sample block can be heated as high as 100 °C. Before removing the RNase P plate or RNase P TLDA, be sure to wait until the sample block reaches room temperature.

**7.** Proceed to "Analyzing the Instrument Verification Run" below.

# Analyzing the Instrument Verification Run

- 1. Click (or select Analysis > Analyze). The SDS software analyzes the run data. A status bar at the bottom of the plate document window indicates progress; the status bar disappears when the analysis is complete.
- **2.** View the results:
  - **a.** In the Plate Grid, click the box above the A label to select all wells.



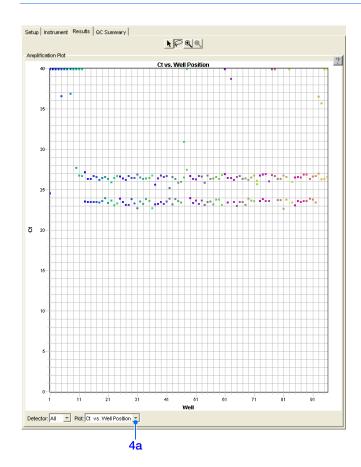
- **b.** Click the **Results** tab to view the plots.
- **3.** Choose from the following:
  - Automatic Ct The SDS software automatically generates baseline values for each well and threshold values for each detector.

Manual Ct – The SDS software calculates baseline and threshold values for a
detector based on the assumption that the data exhibits typical amplification
curves. If you are setting the baseline manually, specify the Start and End
cycles.

**IMPORTANT!** After analysis, you must verify that the baseline and threshold were called correctly for each well by clicking on the well in the Plate Grid and viewing the resulting plots. For more information about manually adjusting the baseline and threshold settings, refer to the *Sequence Detection Systems Software version 2.3 Online Help (SDS Online Help)*.

- **4.** If necessary, remove outliers:
  - **a.** Determine the outlier well positions.

**Note:** The  $C_T$  vs. Well Position view in the Results tab can help you determine the outlier well positions. To access this view, click the **Plot** drop-down list at the bottom of the Amplification Plot.





**b.** In the Plate Grid, select the outlier wells.

**Note:** To select more than one well at a time, hold down the **ctrl** key while selecting the wells.

The number of wells that can be removed to pass specifications is shown in the table below.

Format	Maximum No. of Wells that Can be Removed from Each Set						
	Unknowns	Standards					
Standard 96-well RNase P Plate	6	1					
Fast 96-well RNase P Plate	6	1					
Standard 384-well RNase P Plate	10	2					
RNase P TLDA	4	4					

- c. Click the **Setup** tab, then select the **Omit Wells(s)** check box. A red × appears in each of the selected wells in the Plate Grid.
- **5.** Click (or select **Analysis** > **Analyze**) to reanalyze the data.
- **6.** Proceed to "Verifying Instrument Performance" on page 47.

# **Verifying Instrument Performance**

# Specification Calculation

During analysis, the SDS software generates a standard curve from the averaged threshold cycle ( $C_T$ ) values of the replicate groups of standards, then calculates the concentration of the two unknown populations using the standard curve. To complete the verification, the average copy number (mean quantity) and standard deviation for the unknown populations are entered into the following formula to assess the instrument performance:

$$(Qty Unk_{(high copy)}) - 3 (Qty \sigma_{(high copy)}) > (Qty Unk_{(low copy)}) + 3 (Qty \sigma_{(low copy)})$$

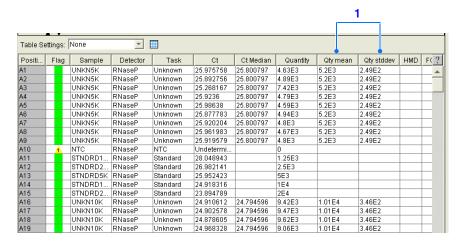
### where:

		Unknown Populations						
Calculation Term	Value	Standard 96- well RNase P Plate	Fast 96-well RNase P Plate	Standard 384- well RNase P Plate	RNase P TLDA			
Qty Unk (high copy)	Average quantity of high copy number	10,000	10,000	10,000	1600			
Qty σ <sub>(high copy)</sub>	Quantity standard deviation of high copy number							
Qty Unk (low copy)	Average quantity of low copy number	5000	5000	5000	800			
Qty σ <sub>(high copy)</sub>	Quantity standard deviation of low copy number							

The calculation states that the average copy number of the larger (10,000 for reaction plates or 1600 for TLDAs) unknown population minus three standard deviations is greater than the average copy number of the smaller (5000 for reaction plates or 800 for TLDAs) unknown population plus three standard deviations. If the calculation is true, then the instrument passes the validation specification, verifying instrument performance.

# Calculating the Results

1. In the Results Table of the plate document window, locate and record the average copy number (Qty mean) and standard deviation (Qty std dev) for the two unknown populations.



**2.** Insert the values into the specification calculation:

$$(Qty\ Unk_{(high\ copy)}) - 3\ (Qty\ \sigma_{(high\ copy)}) > (Qty\ Unk_{(low\ copy)}) + 3\ (Qty\ \sigma_{(low\ copy)})$$

- **3.** Determine whether the specification calculation is true or false.
  - If the calculation is true, the 7900HT Fast System passes the validation specification. Instrument performance is verified.
  - If the calculation is false, the 7900HT Fast System fails the validation specification. The instrument performance cannot be verified.

Notes			

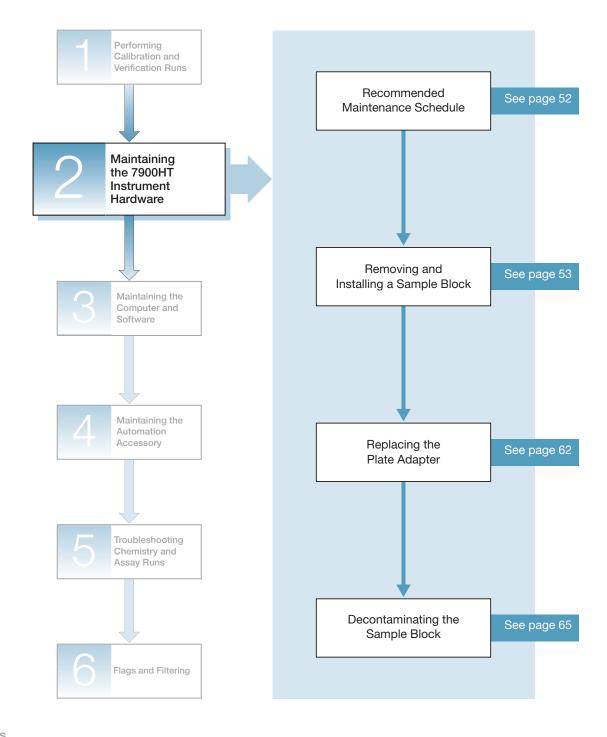
# Chapter 1 Performing Calibration and Verification Runs Performing an Instrument Verification Run

Notes\_\_\_\_\_\_

50 7900HT Fast System Maintenance and Troubleshooting Guide

2

# Maintaining the 7900HT Instrument Hardware



Notes\_\_\_\_

# **Recommended Maintenance Schedule**

To ensure optimal performance of your 7900HT Fast System, perform the following tasks as indicated in the table below.

Schedule	Maintenance Task
Weekly	Cycle the computer and instrument power (power off, then power on the computer and instrument)
Su M T W Th F S	Wipe instrument surfaces with a lint-free cloth
Week (7 Days)	IMPORTANT! Never use organic solvents to clean the 7900HT Fast System.
As needed	Remove and install a sample block (see page 53)
	Replace the plate adapter (see page 62)
	Decontaminate the sample block (see page 65)

Notes			

# Removing and Installing a Sample Block

### When to Perform

You need to remove/install a sample block when you:

- Decontaminate the wells of the sample block (see page 65)
- · Change sample blocks

# Before You Begin

# Time Required

15 minutes. The time will be longer if you have to perform background or pure dye calibrations, or if you have an Automation Accessory.

# Materials Required

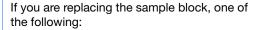
You need the materials listed below to perform this procedure.

5/32 inch hex key

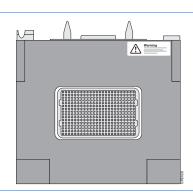
**Note:** Some instruments require the 5/32-inch hex key to remove the sample block locking bar. Other instruments have only a thumbscrew securing the sample block locking bar to the instrument chassis.



**Note:** Some instruments may require a crescent wrench rather than the 5/16-inch hex key.



- 7900HT System Standard 384-Well Block
- 7900HT System Standard 96-Well Block
- 7900HT System Fast 96-Well Block
- 7900HT System TaqMan<sup>®</sup> Low Density Array Block



# Workflow

### To remove/install the sample block:

**1.** Be sure all required upgrades to the SDS software and 7900HT instrument firmware are installed (see page 83).

**IMPORTANT!** Failure to upgrade the software and firmware can make the instrument inoperable or result in damage to instrument components.

- **2.** Obtain the tools and materials required (shown above).
- **3.** Review the sample block handling information (see page 55).



- **4.** Remove the sample block (see page 56).
- **5.** Reinstall the same sample block or install a new sample block (see page 58).
- **6.** After the sample block is successfully installed, perform a background calibration (see page 3) to verify that the sample block:
  - Is connected and working properly
  - Contains no contaminants that will interfere with fluorescent detection

**IMPORTANT!** Perform a background calibration after installing the sample block, even if you are reinstalling the same sample block or replacing it with another block of the same format.

### If you installed a new, uncalibrated sample block, also do the following:

- 7. Change the plate adapter (see page 62).
- **8.** Perform a pure dye calibration to create the pure dye calibration values for the new sample block (see page 17).
- **9.** Perform a verification run to confirm the proper operation of the new sample block (see page 32).

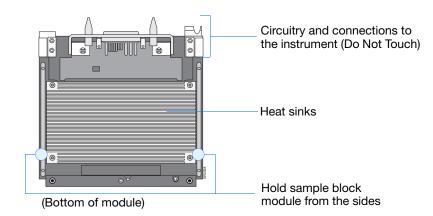
### If you are using an Automation Accessory, also do the following:

- **10.** If you changed sample block formats (for example, replacing a Standard 384-Well Block with a Fast 96-Well Block), adjust the plate sensor switch on the Zymark<sup>®</sup> Twister Microplate Handler arm for the new plate format (see page 92).
- **11.** Align the Plate Handler for the new plate format (see page 97).
- **12.** Align the fixed-position bar code reader for the new plate format (see page 110).

Notes			

# Handling the Sample Block

The interchangeable sample blocks are delicate pieces of equipment containing several fragile components that can break if handled improperly. The figure below shows the correct locations for handling the interchangeable sample blocks.



# Removing the Sample Block

**1.** Confirm the function of the current module:

**Note:** If you do not have the SDS Automation Controller software, skip to step 2 below.

- a. Double-click on the desktop (or select Start > All Programs > Applied Biosystems > SDS 2.3 > SDS Automation Controller 2.3) to start the Automation Controller Software.
- **b.** Select the **Run Status** tab. The module is operating normally if the software is receiving a temperature reading.
- **c.** Click **Open/Close Door** to rotate the instrument tray to the OUT position.
- **d.** Select **File** > **Exit** to close the Automation Controller Software.
- **2.** Power off and unplug the 7900HT instrument.



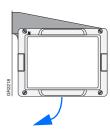
The instrument must be unplugged and powered off at all times during the following procedure. Failure to comply can result in serious physical injury to the user or damage to the instrument.

**3.** Wait 20 to 30 min for the heated cover to cool.

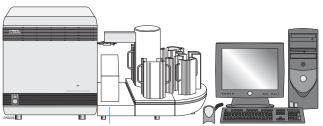


During instrument operation, the temperature of the sample block can be as high as 100 °C. Before performing this procedure, wait until the sample block reaches room temperature.

**4.** If the instrument tray is in the OUT position (outside of the instrument), push it into the instrument to provide an open workspace.



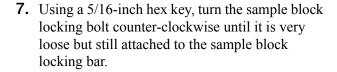
**5.** If using a Plate Handler, remove the covers for the fixed-position bar code reader and the underlying platform.



Fixed-position bar code reader and underlying platform covers

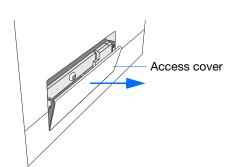
**6.** Push the instrument tray inside the instrument, then remove the thermal cycler access cover to permit access to the sample block.

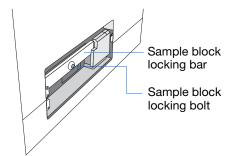
**IMPORTANT!** The thermal cycler access cover is secured to the instrument by non-locking pins and may require force to remove it (no tools are required).

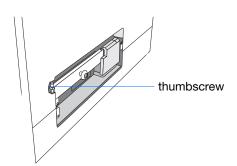


**IMPORTANT!** Some instruments may require the use of an adjustable crescent wrench to loosen the sample block locking bolt.

**8.** Loosen the thumbscrew securing the sample block locking bar to the instrument chassis (may be a 5/32-inch hex bolt on some instruments).

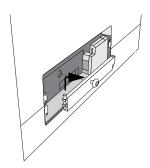






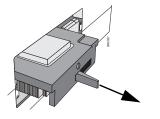


**9.** Lift the sample block locking bar up and out of the instrument.



# **10.** Remove the sample block from the instrument:

- **a.** Rotate the release lever at the base of the sample block 90 degrees.
- b. Being careful not to damage the heat sinks on the bottom of the sample block, slide the sample block out of the instrument and place it on a clean, level surface.

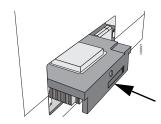


# Installing a Sample Block

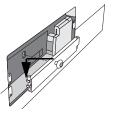
**IMPORTANT!** Before installing the sample block, perform all required upgrades to the SDS software and instrument firmware. Failure to upgrade the software and firmware can render the instrument inoperable or result in damage to instrument components.

- **1.** Load the sample block into the instrument compartment:
  - a. Being careful not to damage the heat sinks on the bottom of the sample block, rest the sample block on the metal runners on either side of the instrument bay.

- b. Carefully slide the sample block into the instrument until the front of the block is flush with the rear of the locking bar.
- **c.** After it is seated, firmly press on the sample block to ensure a good connection.

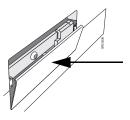


**2.** Replace the sample block locking bar.



- **3.** Tighten the thumbscrew (from step 8 on page 57) to secure the sample block locking bar to the instrument chassis (may be a 5/32-inch hex bolt).
- **4.** Using the 5/16-inch hex key, turn the sample block locking bolt clockwise until it is flush with the locking bar.
- **5.** Again, press on the right and left sides of the front surface of the sample block to ensure that it is seated securely.
- **6.** Replace the thermal cycler access cover:
  - **a.** Fit the lip at the bottom of the access cover over the lower edge of the bay.
  - **b.** Push the cover towards the instrument until it snaps into place.

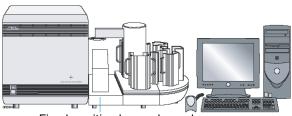
**IMPORTANT!** You must reinstall the thermal cycler access cover before powering on the instrument. Failure to do so prevents power from being applied to the thermal cycler electronics, heated clamp motor, and laser safety interlocks.



# Chapter 2 Maintaining the 7900HT Instrument Hardware Removing and Installing a Sample Block

2

7. If using a Plate Handler, replace the covers for the fixed-position bar code reader and the underlying platform (removed in step 5 on page 57).



Fixed-position bar code reader and underlying platform covers

- **8.** Plug in and power on the 7900HT instrument.
- **9.** Confirm the function of the installed sample block:

**Note:** If you do not have the SDS Automation Controller software, skip this step.

a. Double-click on the desktop (or select Start > All Programs > Applied
 Biosystems > SDS 2.3 > SDS Automation
 Controller 2.3) to start the Automation
 Controller Software.

Notes\_\_\_\_\_

**b.** Select the **Run Status** tab. The sample block is operating normally if the software is receiving a temperature reading.

Does the software display temperatures?	Then
Yes	the installation is successful.
	The presence of temperature readings confirm that the 7900HT instrument successfully established the connection to the new sample block.
No	the 7900HT instrument is unable to establish communication with the new sample block.
	To troubleshoot the problem:
	Power off and unplug the 7900HT instrument.
	2. Remove the thermal cycler access cover.
	<ol><li>Press on the right and left sides of the front plate of the sample block to ensure that it is seated securely.</li></ol>
	4. Reinstall the thermal cycler access cover.
	5. Repeat step 8 and step 9 on page 60 until you hear a high-pitched tone confirming communication between the instrument and sample block.

# Replacing the Plate Adapter

### When to Perform

Replace the 7900HT instrument plate adapter after:

• Changing the sample block format (for example, replacing a Standard 384-Well Block with a Fast 96-Well Block)

**Note:** The sample block must be used with a corresponding plate adapter (that is, both the sample block and the plate adapter must have the same plate format).

# Before You Begin

**Time Required** 

5 to 10 minutes

# Materials Required

You need the materials listed below to perform this procedure.

Flat-blade screwdriver

One of the following:

- 384-Well Plate Adapter
- 96-Well Plate Adapter
- Fast 96-Well Plate Adapter
- TaqMan<sup>®</sup> Low Density Array Adapter

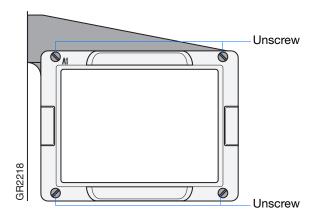
### Workflow T

### To replace the plate adapter:

- **1.** Obtain the tools and materials required (shown above).
- **2.** Replace the plate adapter (see page 63).

# Replacing the Plate Adapter

- **1.** If the instrument tray is inside the 7900HT instrument, move the instrument tray to the OUT position:
  - a. Double-click on the desktop (or select Start > All Programs > Applied Biosystems > SDS 2.3 > SDS 2.3) to start the SDS software.
  - **b.** Click  $\Box$  (or select **File** > **New**).
  - c. In the New Document dialog box, click | ok |.
  - **d.** In the new plate document, select the **Instrument > Real-Time** tabs.
  - e. Click Open/Close
  - **f.** The instrument tray rotates to the OUT position.
  - g. Select **File** > **Exit** to close SDS software.
- **2.** Remove the four screws attaching the plate holder to the plate arm.



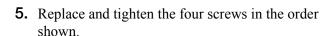
**3.** Remove the plate adapter from the instrument tray.

**Note:** If changing sample block formats (for example, replacing a Standard 384-Well Block with a Fast 96-Well Block), store the plate adapter with the sample block of the same format.

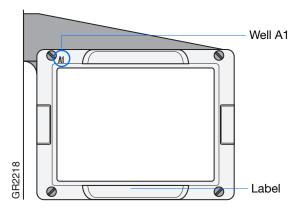


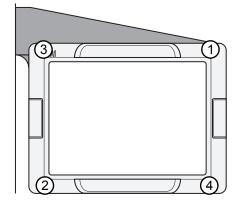
**4.** Place the new plate adapter into the instrument tray with the A1 label in the rear-left corner.

**IMPORTANT!** Make sure to install the correct version of the plate adapter for the plate format you intend to use. The plate adapters are labeled with the format they support.



**IMPORTANT!** The order in which the screws are tightened is important to ensure proper alignment of the plate to the sample block inside the 7900HT instrument.





# **Decontaminating the Sample Block**

# Overview

This section describes how to decontaminate the wells of a sample block. The procedure eliminates residual PCR-related products, including fluorescent labeled TaqMan® probes.

# Using Other Methods

**IMPORTANT!** If you plan to use a decontamination method other than the one in this guide, check with Applied Biosystems first to ensure that the method will not damage the sample block or the 7900HT instrument.

### When to Perform

Decontaminate the sample block as often as needed.

**Note:** Decontamination is generally performed to resolve problematic background calibrations, where one or more wells consistently exhibit abnormally high signals indicating the presence of a fluorescent contaminant.

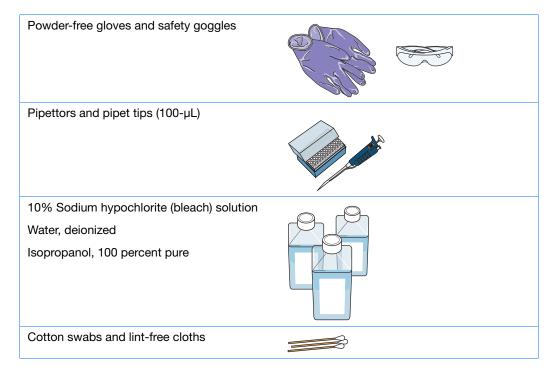
# Before You Begin

Time Required

30 minutes. The time may vary, depending on the extent of the contamination.

# Materials Required

You need the materials listed below to perform this procedure.





# Workflow To decontaminate the sample block:

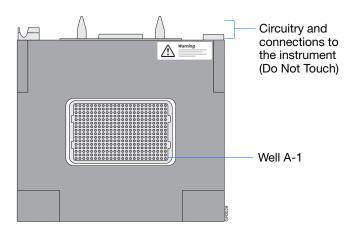
- **1.** Remove the sample block from the 7900HT instrument (see page 53).
- **2.** Obtain the tools and materials required for decontamination (shown above).
- **3.** Decontaminate the sample block:
  - For a Standard 384-Well Block, Standard 96-Well Block, or Fast 96-Well Block, see page 67
  - For a TaqMan® Low Density Array Block, see page 70
- **4.** Replace the sample block (see page 58).
- **5.** Perform a background calibration to confirm that the contamination has been removed (see page 3).

Notes			

# Decontaminating a Sample Block

This procedure applies to a:

- Standard 384-Well Block
- Standard 96-Well Block
- Fast 96-Well Block
- 1. Suspected contaminated wells are those that give a signal >4000 FSU during a background calibration (see page 14). Using the SDS software, isolate the suspected contaminated wells as explained on page 15.
- **2.** Locate the suspected contaminated wells on the sample block, using the figure at right as a guide.





- **3.** Rinse (pipet and remove) each contaminated well with three treatments of deionized water at the appropriate volume for the sample block:
  - 40 μL per well for a Standard 384-Well Block
  - 150 μL per well for a Standard 96-Well Block
  - 55 μL per well for a Fast 96-Well Block

**Note:** Absolute isopropanol can be substituted for water in the third treatment.

# WARNING CHEMICAL HAZARD.

**Isopropanol** is a flammable liquid and vapor. It may cause eye, skin, and upper respiratory tract irritation. Prolonged or repeated contact may dry skin and cause irritation. It may cause central nervous system effects such as drowsiness, dizziness, and headache, etc. Please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

**4.** Using a cotton swab, scrub inside of each contaminated well.

**Note:** If you are decontaminating a Standard 384-Well Block, remove some of the cotton from the swab tip; this allows the swab to reach the bottom of the wells.

**5.** Using a lint-free cloth, absorb the excess deionized water or isopropanol from the wells.

Notes			

- **6.** Pipet the appropriate volume of 10% bleach solution into each suspected contaminated well:
  - 40 μL per well for a Standard 384-Well Block
  - 150 µL per well for a Standard 96-Well Block
  - 55 μL per well for a Fast 96-Well Block

WARNING CHEMICAL HAZARD.

Sodium hypochlorite (bleach) is a liquid disinfectant that can be corrosive to the skin and can cause skin depigmentation. Please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

- **7.** Allow the sample block to sit for 3 to 5 min.
- **8.** Using a pipet, remove the bleach solution from the wells.
- **9.** Rinse (pipet and remove) the wells with deionized water at the appropriate volume for the sample block:
  - 40  $\mu L$  per well for a Standard 384-Well Block
  - 150 µL per well for a Standard 96-Well Block
  - 55 μL per well for a Fast 96-Well Block
- **10.** Using a lint-free cloth, absorb the excess deionized water from the wells.

- N	ь	$\overline{}$	4.	~
1	A (	U	U	E:

# Decontaminating a TLDA Block

This procedure applies to the TaqMan Low Density Array Block.

**1.** Using a lint-free cloth, wipe the surface of the sample block (grey aluminum) with 10% bleach.

WARNING CHEMICAL HAZARD.
Sodium hypochlorite (bleach) is a liquid disinfectant that can be corrosive to the skin and can cause skin depigmentation. Please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

- **2.** Wipe the sample block three times with distilled water.
- **3.** Wipe the sample block with isopropanol.

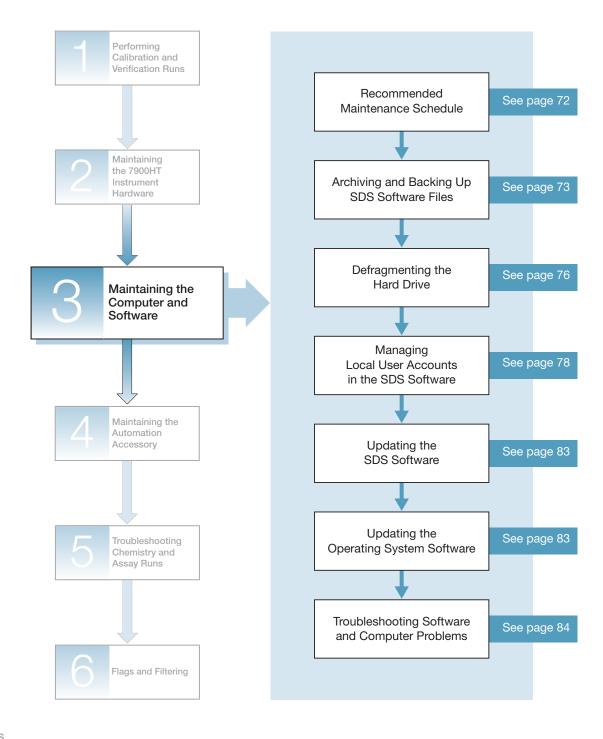
# WARNING CHEMICAL HAZARD. Isopropanol is a flammable liquid and vapor. It may cause eye, skin, and upper respiratory tract irritation. Prolonged or repeated contact may dry skin and cause irritation. It may cause central nervous system effects such as drowsiness, dizziness, and headache, etc. Please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

**4.** Allow the sample block to air dry.

Notes		

# 3

# Maintaining the Computer and Software



# **Recommended Maintenance Schedule**

To ensure optimal performance of your 7900HT Fast System, perform the following tasks as indicated in the table below.

Schedule	Maintenance Task
Weekly	Archive or backup SDS plate document files (see page 73)
Su M T W Th F S  Week (7 Days)	
Monthly  Su M T W Th F S  Month (30 Days)	Defragment the computer hard drive (see page 76)
As needed	<ul> <li>Manage local user accounts in the SDS software (see page 78)</li> <li>Update the SDS software (see page 83)</li> <li>Update the operating system software (see page 83)</li> </ul>

# Archiving and Backing Up SDS Software Files

### When to Perform

Weekly

# Archiving SDS Software Files

To conserve space on the computer hard drive, SDS software files can be archived using a data compression utility. A compression utility typically archives files by encoding the binary data they contain algorithmically, thereby reducing the size of a file.

Several commercially available compression utilities are available. PKZIP and \*.arc are archive formats common to the Microsoft Windows® operating system.

# Backing Up SDS Software Files

Applied Biosystems strongly recommends that you back up the data generated by your 7900HT Fast System for two reasons:

- Backing up data protects against potential loss of data caused by an unforeseen failure of the computer or its hard drive(s).
- Backing up older data and removing it from the computer hard drive conserves space on the hard drive and optimizes performance.

### Choosing a Backup Storage Device

Applied Biosystems recommends the use of one or more backup storage devices to prevent potential loss of data caused by unforeseen failures of the computer or its hard drive(s).

The CD-RW drive of the computer can serve as the backup storage device for your system. By saving your \*.sds and \*.sdt files to one or more writable CDs on a weekly basis, you can effectively backup the data generated by your 7900HT Fast System.

# Developing a Data Management Strategy

Applied Biosystems recommends developing a strategy for dealing with the files produced by the SDS software. During a single day of real-time operation, the 7900HT System can generate over 200 MB of data. Without a strategy for distributing and archiving SDS software files, the 7900HT instrument can easily fill the hard drive of the computer within just a few weeks of operation.

If you are performing real-time experiments on your 7900HT Fast System, check the amount of available space on your hard drive weekly. When the hard drive is within 20% of it's capacity, transfer the older data to a backup storage device.

**Note:** Real-time experiments include standard curve (AQ) assays and comparative C<sub>T</sub> (RQ) assays, as well as any amplification (AQ) runs you might perform on the 7900HT System for allelic discrimination assays or plus/minus assays.

To successfully manage the information produced by the 7900HT System, you should have a basic understanding of how data is collected, stored, and processed throughout the operation of the instrument. See "Modes of Operation" on page 74.

Notes		

# **Modes of Operation**

# System Operation and Dataflow

Data management strategy is a crucial element of successfully integrating the Applied Biosystems 7900HT Fast Real-Time PCR System into a laboratory workflow. During a single 24-hour period of real-time operation, the 7900HT instrument can produce more than 200 MB of data. To successfully manage the information produced by the 7900HT System, you should have a basic understanding of how data is collected, stored, and processed throughout the operation of the instrument.

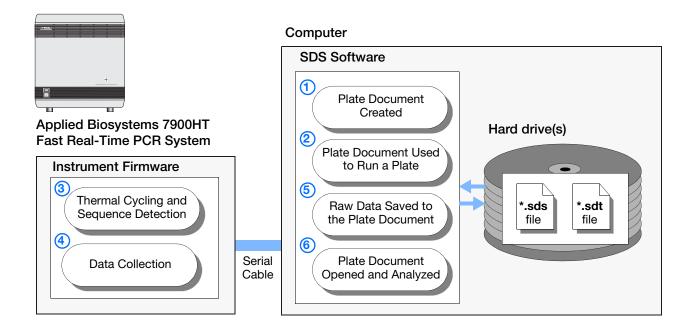
# Standard Modes of Operation

The 7900HT Fast System has two standard modes of operation, depending on the accessories purchased with the base instrument:

- Stand-Alone Operation (see below)
- Automated Operation (see page 75)

### **Stand-Alone Operation**

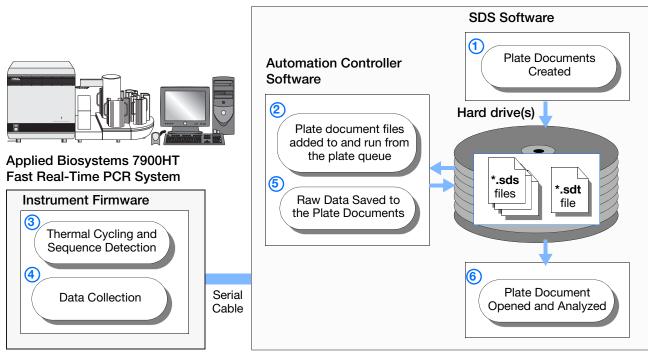
In this mode (the most basic configuration of the 7900HT Fast System), the 7900HT instrument is used without any additional components. In this configuration (see the figure below), a technician runs plates individually using the SDS software. The computer stores all data for instrument operation and provides the software for analyzing the run data.



### **Automated Operation**

In this mode, the 7900HT Fast System uses an Automation Accessory (Zymark® Twister Microplate Handler and fixed-position bar code reader) to provide high-throughput operation suitable for small- to medium-scale studies. With this configuration (see the figure below), a technician can use the Automation Controller Software to run batches of plates unattended. As in Stand-Alone mode, the computer stores all data for instrument operation and provides the software for analyzing the run data.

### Computer



# **Defragmenting the Hard Drive**

### When to Perform

- At least once every month
- Before fragmentation reaches 10% (or when warned by the Windows operating system)

# Purpose: Why Defragment the Hard Drive?

As the 7900HT Fast System is used and files are deleted and created, the free space on the computer hard drive eventually is split into increasingly smaller blocks. Consequently, as the SDS software creates new files and extends old ones, the computer cannot store each file in a single block. Instead, the system will 'fragment' the files by scattering their component pieces across different sectors of the hard drive.

The fragmentation of SDS files decreases the performance of both the SDS software and the computer operating system. As the hard drive becomes fragmented, programs take greater time to access files because they must perform multiple seek operations to access the fragments. Defragmentation utilities defragment broken files by combining their component pieces at a single location on the hard drive, thereby optimizing system performance.

Notes		

# Defragmenting the Hard Drive

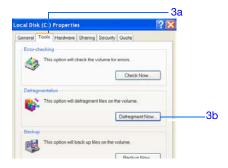
- 1. In the Windows desktop, select **y** start **My Computer**.
- **2.** In the My Computer window, right-click a hard drive and select **Properties**.

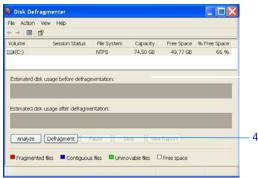
- **3.** In the Local Disk Properties dialog box:
  - a. Select the Tools tab.
  - b. Click DefragmentNow... to open the Disk Defragmenter dialog box.
- 4. Click Defragment.



- **6.** In the Local Disk Properties dialog box, click ox .
- **7.** Repeat steps 2 through 6 for the remaining drives on the computer.







# Managing Local User Accounts in the SDS Software

Managing local user accounts includes the following:

- Enabling or disabling the SDS software login function (page 79)
- Adding user accounts (page 79)
- Editing user accounts (page 81)
- Deleting user accounts (page 82)

### Non-Enterprise Mode Only

This section applies to the SDS software in non-Enterprise mode (that is, the SDS software does not include the SDS Enterprise Database). For versions of the SDS software that do include the SDS Enterprise Database, user accounts are managed with the SDS User Account Manager software. For more information, see the SDS Enterprise Database for the Applied Biosystems 7900HT Fast Real-Time PCR System Administrators Guide.

When to Perform As needed

Notes		

# **Enable/Disable Login**

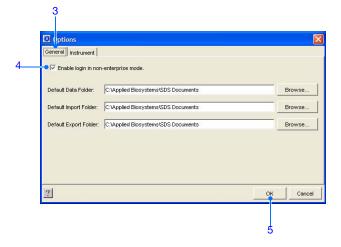
Enabling the login function allows the SDS software to capture user information so that the system can be customized for each user (user preferences).

When the login is enabled, the Login dialog box appears when the SDS software is started.



### To enable or disable the login:

- Double-click on the desktop (or select Start > All Programs > Applied Biosystems > SDS 2.3 > SDS 2.3) to start the SDS software.
- **2.** Select **Tools ▶ Options** to open the Options dialog box.
- **3.** Click the **General** tab.
- **4.** Select or deselect the **Enable login in non-enterprise mode** check box, as desired.
- **5.** Click **OK** to save the changes and close the dialog box.

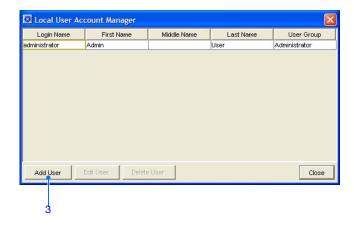


### Add a User Account

Double-click on the desktop (or select Start > All Programs > Applied Biosystems > SDS
 2.3 > SDS
 2.3) to start the SDS software.



- 2. Select Tools > Local User Account Manager to open the Local User Account Manager dialog box.
- 3. Click Add User.



- **4.** Complete the Add Local User dialog box:
  - a. User Name Enter a user name. The user name must begin with a letter of the alphabet and cannot exceed 30 characters.
  - b. Password/Confirm Password Enter a password, then re-enter it to confirm. The password must begin with a letter of the alphabet and must be between 6 and 30 characters long.

Note: Users can change their own passwords at a later time by selecting Tools > Change Password.

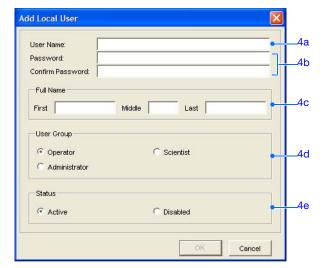
- **c.** Full Name pane Enter the user's First, Middle, and Last names. The names cannot exceed 50 characters.
- d. User Group pane Select an appropriate group for the user: Operator,
   Administrator, or Scientist.

**Note:** This information is *not* used for access control purposes.

**e.** Status pane – Select an appropriate status for the user: **Active** or **Disabled**.

**Note:** If a user has the Disabled status, the user cannot log into the software.

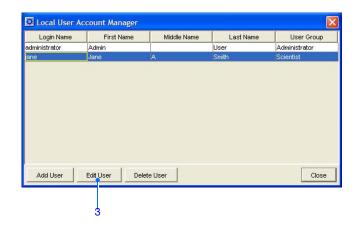
**5.** Click **OK** to save the changes and close the dialog box.



3

### Edit a User Account

- Double-click on the desktop (or select Start > All Programs > Applied Biosystems > SDS
   2.3 > SDS
   2.3) to start the SDS software.
- 2. Select Tools > Local User Account Manager to open the Local User Account Manager dialog box.
- **3.** Select the user account you want to edit, then click **Edit User**.



- **4.** Complete the Edit Local User dialog box:
  - a. Full Name pane Enter the user's First, Middle, and Last names. The names cannot exceed 50 characters.
  - b. User Group pane Select an appropriate group for the user: Operator,Administrator, or Scientist.

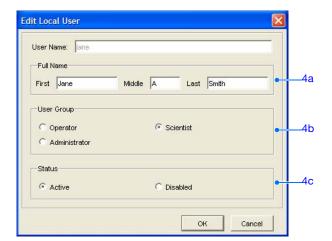
**Note:** This information is *not* used for access control purposes.

**c.** Status pane – Select an appropriate status for the user: **Active** or **Disabled**.

**Note:** If a user has the Disabled status, the user cannot log into the software.

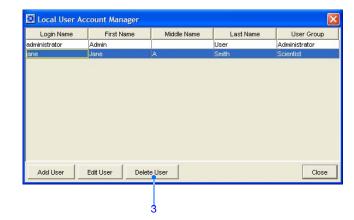
Note: You cannot edit a user's name or password. Users can change their own passwords at a later time by selecting Tools ▶ Change Password.

**5.** Click **OK** to save the changes and close the dialog box.



### Delete a User Account

- Double-click on the desktop (or select Start > All Programs > Applied Biosystems > SDS 2.3 > SDS 2.3) to start the SDS software.
- 2. Select Tools ➤ Local User Account Manager to open the Local User Account Manager dialog box.
- **3.** Select the user account you want to delete, then click **Delete User**.



**4.** At the prompt, click **Yes**. The user account is removed from the Local User Account Manager.



**5.** Click **Close** to close the dialog box.

# **Updating the SDS Software**

### When to Perform

As needed

# Upgrading the SDS Software

Applied Biosystems continually develops the SDS software to provide increased functionality and reliability of the 7900HT Fast System. As updates become available, Applied Biosystems sends notifications of the upgrades to all Applied Biosystems 7900HT Fast System customers.

Review all documentation accompanying the software upgrade (such as installation notes or user bulletin). The updated version of the software may contain new features that require special consideration.

**Note:** Applied Biosystems service engineers perform regular SDS software updates during planned maintenance visits.

# Reinstalling the SDS Software

On rare occasions, when a piece of the SDS software becomes corrupt, it may be necessary to reinstall the software. In the event that the software must be reinstalled, observe the following guidelines:

- Unless instructed to do otherwise, remove the SDS software using the uninstall utility. Do not delete the program folder from the Program Files directory.
- Backup all data before reinstalling the SDS software.
- Reinstall the SDS software under a user login that has administrator privileges on the computer.
- Unless instructed to do otherwise, reinstall the SDS software to the same directory as the previous installation.

# **Updating the Operating System Software**

When to Perform Windows Service Pack Updates Do not upgrade the operating system of the computer connected to the 7900HT Fast System unless instructed to do so by an Applied Biosystems representative. New versions of the Windows operating system can conflict with the SDS software and render the instrument inoperable.

If you want to install a service pack to update the operating system, check the release notes that are installed with the SDS software for compatibility issues.

**Note:** Applied Biosystems service engineers maintain the operating system software as part of planned maintenance visits. During the visit, an engineer will update the computer operating system as upgrades become available and are validated by Applied Biosystems.

Notes		

# **Troubleshooting Software and Computer Problems**

# Troubleshooting Table

Table 3-1 Troubleshooting software and computer problems

Observation	Possible Cause	Recommended Action
SDS software will not start The software crashes/freezes	Incorrect start-up sequence	Follow the solutions listed until the symptom goes away.
the computer or displays an error message	<ul><li>Corrupted software</li><li>Computer hardware failure</li></ul>	1
	Operating System (OS) corruption	Power off the 7900HT instrument.     Check cable connections.
	Loose bar code reader cable	Restart the computer and log on to the computer.
		Power on the 7900HT instrument.
		5. Start the SDS software.
		2
		Restart the computer and log on to your computer.
		7. Reinstall the SDS software.
		8. Start the SDS software.
		3
		Contact Applied Biosystems Service for OS problems or if the computer will not boot up at all. You may have to reload the OS from the CDs.
		4
		Contact Dell for troubleshooting the computer hardware.
Communication error	Cables are connected incorrectly	Check cable connections and COM port setup. See Appendix B on page 153.
Thermal cycler errors	Sample block module not fully engaged	Reseat the sample block module. See "Removing and Installing a Sample Block" on page 53.
Automation Controller Software cannot find a plate document file	File not in correct location	Remove file entry from plate queue and add the file to the plate queue again.

Table 3-1 Troubleshooting software and computer problems

Observation	Possible Cause	Recommended Action
Dialog box does not respond to mouse clicks or key strokes	Java Runtime Error	Click the close box of the dialog box to close it.
Run will not start	No calibration file	Perform background and pure dye calibrations.
	No background data in calibration file (background calibration has not been performed)	See "Performing a Background Calibration" on page 3 and "Performing a Pure Dye Calibration" on page 17.
	No pure dye data in calibration file (pure dye run has not been performed)	
	Calibration file does not contain pure dye data for a dye used on the plate document	
	Calibration file was created on another instrument	
	Disk drive containing the plate document has less than 50 MB of free space	Check the capacity of the destination drive. If less than 50 MB of free space remains, remove or archive existing data files. See "Archiving and Backing Up SDS Software Files" on page 73.
	Heated cover cannot reach running temperature because no plate loaded	Open the instrument tray and check that the instrument contains a plate.
	Instrument tray contains a plate	
	Output stack contains a plate or plates	Remove all plates from the output stack of the Plate Handler before starting the queue.
Computer is slow when analyzing data, opening or closing dialog boxes, and other software processes	Hard drive is fragmented	Defragment the hard drive. See "Defragmenting the Hard Drive" on page 76.
	Hard drive is almost full	Remove or archive existing data files. See "Archiving and Backing Up SDS Software Files" on page 73.



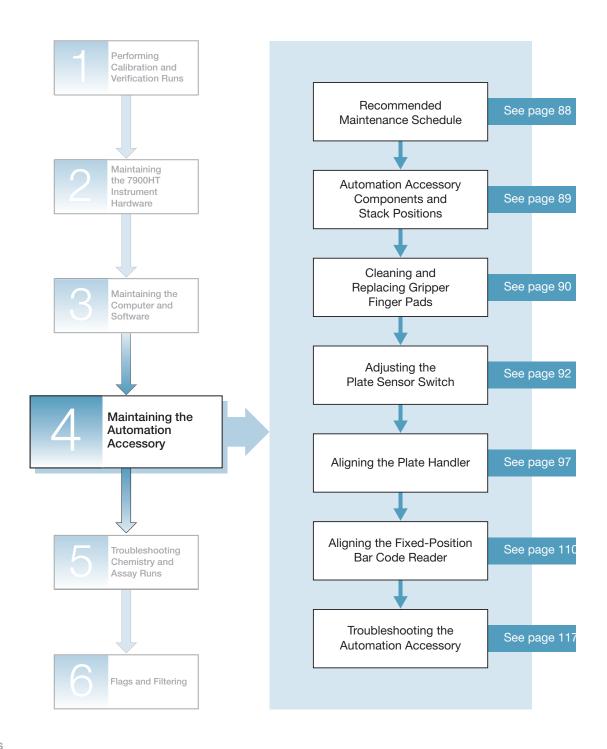
Table 3-1 Troubleshooting software and computer problems

Observation	Possible Cause	Recommended Action
The computer will not logon to the Windows operating system	Logon window does not appear	Restart the computer and log on to your computer.
System:	You are not logged on as the <b>Administrator</b>	<ol> <li>Log off of your computer.</li> <li>Log on again as the Administrator.</li> </ol>
	After the above solutions have been tried, the problem is still not fixed	Contact Dell for troubleshooting the computer hardware or OS.
The computer will not boot up at all	Cables are not connected or are not	Check the cables.
up at all	seated properly	The boot disk is corrupted.
		Boot directly off of the     Windows operating system     installation CD.
		Boot off of the emergency disk.
		Reload the Windows operating system from the CD.
	After the above solution has been tried, the problem is still not fixed	Contact Dell for troubleshooting the computer hardware.

Notes\_\_\_\_



# Maintaining the Automation Accessory



# **Recommended Maintenance Schedule**

To ensure optimal performance of the Automation Accessory, perform the following tasks as indicated in the table below.

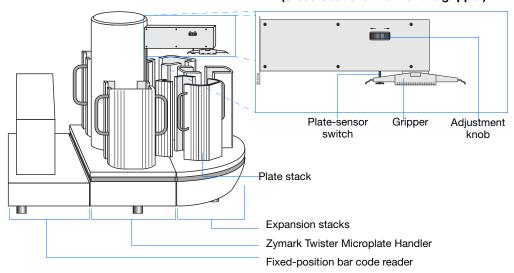
Schedule	Maintenance Task
Monthly	Inspect the gripper finger pads (see page 90)
Su M T W Th F S  Month (30 Days)	
As needed	Clean or replace the gripper finger pads (see page 90)
	<ul> <li>Adjust the plate sensor switch (see page 92)</li> </ul>
	Align the Plate Handler (see page 97)
	Align the fixed-position bar code reader (see page 110)

# **Automation Accessory Components and Stack Positions**

Automation Accessory Components The Automation Accessory includes the Zymark® Twister Microplate Handler and the fixed-position bar code reader.

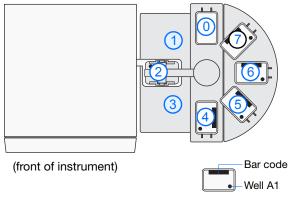
Refer to the figure below for the components discussed in this section.

### (cross-sectional view of the gripper)



### Plate Stack Positions

The Zymark Twister Microplate Handler alignment is performed using the Zymark® Twister Software. The software refers to the positions of the plate stacks differently than the Automation Controller Software. The figure below lists the positions defined by the Zymark Twister Software and the Automation Controller Software equivalents.



Zymark Twister Software	Automation Controller Software
Position 0	Output
Position 1	(unused)
Position 2	Instrument
Position 3	(unused)
Position 4	Stack 1

# Cleaning and Replacing Gripper Finger Pads

### When to Perform

The adhesive used to affix bar code labels to certain brands of microplates can build up on the gripper pads of the Zymark Twister Microplate Handler. Over time, the residue can cause the gripper pads to stick to the microplates while handling them, causing misfeeds.

To prevent buildup, Applied Biosystems recommends that you:

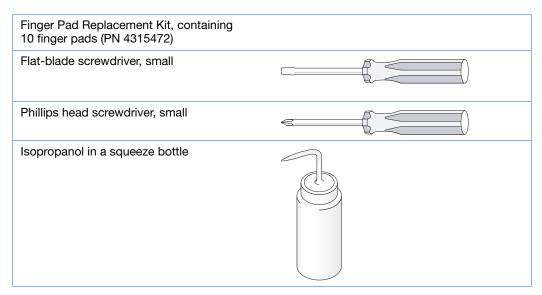
- Inspect the gripper pads monthly
- · Clean or replace the pads as needed

# Before You Begin

Time Required 10 minutes

# Materials Required

You need the materials listed below to perform this procedure.



# Workflow To clean and replace the gripper finger pads:

- 1. Obtain the tools and materials required (shown above).
- **2.** Clean (see below) or replace (see page 91) the gripper finger pads.

4

# Cleaning the Finger Pads

1. Wipe each pad thoroughly with Isopropanol until the residue has been removed.

WARNING CHEMICAL HAZARD. Isopropanol is a flammable liquid and vapor. It may cause eye, skin, and upper respiratory tract irritation. Prolonged or repeated contact may dry skin and cause irritation. It may cause central nervous system effects such as drowsiness, dizziness, and headache, etc. Please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

**2.** If the pads appear rough or the adhesive cannot be removed, replace the pads as described on page 91.

# Replacing the Finger Pads

1. Using a small Phillips-head screwdriver, remove the two small Phillips-head screws from the fingers on each side of the gripper, then remove the fingers.

**Note:** Move the Plate Handler arm into any position where it is easy to access the screws.

**2.** Using a small flat-blade screwdriver, pry the worn finger pads off the fingers.

**Note:** The manufacturer recommends replacing all finger pads at the same time.

**3.** Clean any residual adhesive off the fingers using isopropanol.

WARNING CHEMICAL HAZARD. Isopropanol is a flammable liquid and vapor. It may cause eye, skin, and upper respiratory tract irritation. Prolonged or repeated contact may dry skin and cause irritation. It may cause central nervous system effects such as drowsiness, dizziness, and headache, etc. Please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

- **4.** Remove a replacement finger pad from the paper backing, then place the finger pad on the appropriate finger position.
- **5.** Repeat step 4 for the remaining finger pads.
- **6.** Install the fingers with the fingers pointing down and the finger pads toward the center of the gripper.
- **7.** Insert the screws into the fingers and tighten.

**Note:** The screws do not automatically align the grippers. Make sure that the finger pads are making good contact with the plate when the Plate Handler arm grips a plate.

# Adjusting the Plate Sensor Switch

### When to Perform

Adjust the plate sensor switch (located underneath the arm of the Plate Handler):

• After changing the sample block format (for example, replacing a Standard 384-Well Block with a Standard 96-Well Block)

**Note:** The Automation Accessory should not be used for running Fast 96-well reaction plates.

**Note:** The various plate types have different dimensions, which can affect the way the Plate Handler grips the plates. Adjusting the plate sensor switch after changing sample block formats ensures smooth operation of the Automation Accessory.

• If the Plate Handler is having difficulty sensing plates

# **Before You Begin**

### Time Required

20 minutes

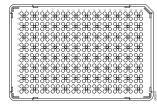
# Materials Required

You need the materials listed below to perform this procedure.

Appropriate plate type for your current sample block; one of the following:

- Standard 384-well reaction plate
- Standard 96-well reaction plate
- TaqMan<sup>®</sup> Low Density Array

**Note:** The Automation Accessory should not be used for running Fast 96-well reaction plates.



Pictured: Standard 384-well reaction plate

### Workflow

To adjust the plate sensor switch:

- 1. Obtain the tools and materials required (shown above).
- **2.** Adjust the plate sensor switch (see page 93).
- **3.** Test the adjustment (see page 95).
- **4.** Confirm the Zymark Twister Software is closed (see page 96).

010	

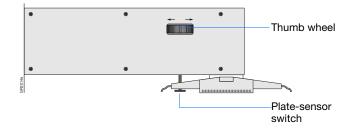
# Adjusting the Plate Sensor Switch

**1.** Power off the Zymark Twister Microplate Handler.



The Zymark Twister Microplate Handler must be powered off at all times during the following procedure. Failure to comply can result in physical injury to the user or damage to the Plate Handler.

**2.** Clear the switch position by turning the thumb wheel all the way to the Up extreme (as indicated on the side panel).



- **3.** Begin the adjustment of the sensor switch:
  - a. Grasp a plate by the sides, making sure not to place pressure in the center of the plate to deform it.
  - **b.** Place the plate between the fingers of the gripper assembly and align it to the middle of the centering device.
  - **c.** While holding the plate in position, slowly turn the thumb wheel to lower the switch onto the plate until the switch:
    - · Contacts the top of the plate, and
    - Emits a soft, audible "clicking" noise

**IMPORTANT!** The sound emitted by the sensor switch is very faint and may be difficult to hear. To make the adjustment easier, place your ear close to the sensor switch while making the adjustment and listen for the switch to engage.



**4.** Remove the plate and listen for the plate sensor switch to disengage.

Did you hear the switch disengage?	Then
No	Move the switch Down a few steps by turning the thumb wheel in the direction indicated on the arm.
	<ul> <li>Replace the plate inside the gripper and listen for the switch to engage:</li> </ul>
	<ul> <li>If you do not hear the switch engage, then remove the plate and repeat steps a and b above.</li> </ul>
	<ul> <li>If you hear the switch engage, remove the plate and continue to step a below.</li> </ul>
Yes	Move the switch Up by turning the thumb wheel one step in the direction indicated on the arm.
	b. Replace the plate and listen for the switch to engage:
	<ul> <li>If you hear the switch engage, remove the plate and repeat steps a and b above.</li> </ul>
	<ul> <li>If you do not hear the switch engage, then you have successfully identified the zero point of the plate sensor switch.</li> </ul>
	<b>Note:</b> At the zero point, one step of the thumb wheel in the Down direction causes the switch to engage.

**5.** After the zero point is established, carefully turn the thumb wheel in the Down direction the number of steps appropriate for your plate type, as indicated below:

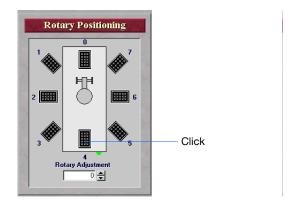
Plate	Turn the thumb wheel in the Down direction
Standard 384-well reaction plate	15 steps
Standard 96-well reaction plate	20 steps
TaqMan Low Density Array	15 steps

**Note:** If you lose count, begin again from step 4 and identify the zero point for the switch.

B. I	_		_	
IVI	$\cap$	١T	$^{\circ}$	c

# **Testing the Adjustment**

- **1.** Place the plate in the input stack 1 of the Plate Handler.
- **2.** Power on the 7900HT Fast System, the Plate Handler, and the computer.
- **3.** Select **Start > Programs > Zymark Twister Plate Handler > Twister** to start the Zymark
  Twister Software.
- **4.** Click **Manual Control** to open the Manual Control dialog box.
- **5.** Click stack **4**. The Plate Handler arm moves over the input stack.



### 6. Click & Find Plate

- If the adjustment was successful, the Plate Handler arm will lower upon the plate until the plate sensor switch engages, confirming the presence of the plate.
- If the Plate Handler arm emits a grinding sound, adjust the plate sensor switch:
- a. In the Zymark Twister Software, click

  to vertical Home to raise the Plate Handler arm.
- **b.** Turn the thumbscrew in the Down direction 10 steps.
- **c.** Repeat step 6 until the Plate Handler arm successfully detects the plate.



- 7. Click Close Gripper, then click Vertical Home
  - If the adjustment was successful, the Plate Handler arm will grasp the plate and remove it from the plate stack.
  - If the Plate Handler arm stops before the gripper fingers are able to contact the plate and fails to grasp or pick up the plate, adjust the plate sensor switch:
  - **a.** Turn the thumbscrew in the Up direction 10 steps.
  - b. Grasp the plate with one hand and, from the Zymark Twister Software, click

    or Open Gripper to release the plate.
  - **c.** Replace the reaction plate into input stack 1 of the Plate Handler.
- **8.** Repeat steps 6 through 7 until the Plate Handler arm successfully retrieves the plate.
- 9. Grasp the plate with one hand and, from the Zymark Twister Software, click Gripper to release the plate.
- **10.** Exit the Zymark Twister Software:
  - a. Click → Main Menu
  - **b.** Click **Quit Application** to close the software.

### Confirm the Zymark Twister Software Is Closed

A defect in the Zymark Twister Software can cause portions of the program to persist in memory even after the software has been closed. Because the Zymark Twister Software conflicts with the SDS software, the residual elements of the Zymark Twister Software must be closed inside the Microsoft Windows® operating system Task Manager before continuing.

- **1.** Press the **Crtl** + **Alt** + **Del** keys in unison to open the Windows operating system Security dialog box.
- **2.** Click **Task Manager** to open the Windows operating system Task Manager dialog box.
- **3.** Click the **Applications** tab.
- **4.** Confirm that the software has closed by looking for the Zymark Twister Software entry in the **Task** list.
  - If the software is completely closed, there is no entry for the Zymark Twister Software.
  - If the software is still running, click the Zymark Twister Software entry to select it, then click **End Task**.
- **5.** Close the Task Manager dialog box.

Notes		

### When to Perform

Align the Zymark Twister Microplate Handler if:

- You move the Applied Biosystems 7900HT Fast Real-Time PCR System
- The Plate Handler is misaligned
   Symptoms that the Plate Handler is out of alignment include:
  - Excessive downward movement of the Plate Handler arm (the arm grinds when grasping or releasing plates)
  - The Plate Handler arm collides with the plate stacks
  - The Plate Handler arm releases plates above the bottom of the plate stacks
  - Plates tip or tilt when placed into the instrument tray by the Plate Handler arm

# Before You Begin

**Time Required** 

30 minutes

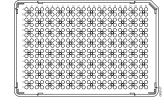
# Materials Required

You need the materials listed below to perform this procedure.

Appropriate plate type for your current sample block; one of the following:

- Standard 384-well reaction plate
- Standard 96-well reaction plate
- TaqMan<sup>®</sup> Low Density Array

**Note:** The Automation Accessory should not be used for running Fast 96-well reaction plates.



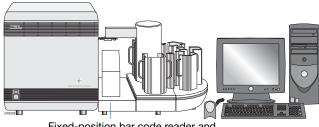
Pictured: Standard 384-well reaction plate

# Workflow To align the plate handler:

- **1.** Obtain the tools and materials required (shown above).
- **2.** Prepare the 7900HT instrument (see page 99).
- **3.** Align input stack 1 (see page 100).
- **4.** Align the Plate Handler to the 7900HT instrument (see page 102).
- **5.** Recheck input stack 1 (see page 104).
- **6.** Define the bottom of the stack (see page 105).
- 7. Define the positions of the remaining stacks (see page 107).
- **8.** Confirm the Zymark Twister Software is closed (see page 108).

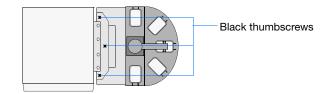
# Preparing the Instrument for the Alignment

**1.** Remove the covers for the fixed-position bar code reader and the underlying platform.



Fixed-position bar code reader and underlying platform covers

**2.** Loosen the three black thumbscrews on the platform connecting the 7900HT Fast System and the Plate Handler base.



- **3.** Power on the 7900HT Fast System, the Plate Handler, and the computer.
- **4.** Move the instrument tray to the OUT position:
  - a. Double-click on the desktop (or select Start > All Programs > Applied Biosystems > SDS 2.3 > SDS Automation Controller 2.3) to start the Automation Controller Software.

Note: If an error dialog box appears reading, Machine calibration values are not valid. Please refer to documentation for calibration process, click \_\_\_ok\_\_|.

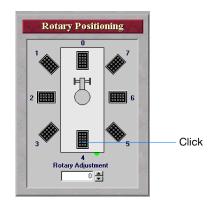
- b. Click Open/Close Door. The 7900HT Fast System moves the instrument tray to the OUT position.
- **c.** Select **File** > **Exit** to close the Automation Controller Software.
- 5. Select Start > Programs > Zymark Twister Plate Handler > Twister to start the Zymark Twister Software.
- **6.** Click **Manual Control** to open the Manual Control dialog box.



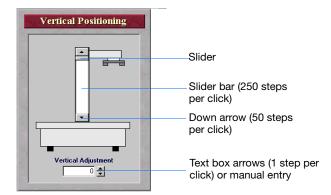
# Aligning Input Stack 1 (Zymark Position 4)

The alignment of input stack 1 (position 4 in the Zymark Twister Software) is the first step in the alignment procedure. This alignment provides the basis for aligning all subsequent stacks on the Plate Handler.

- **1.** Place an empty plate into input stack 1 (Zymark position 4).
- **2.** In the Zymark Twister Software, click position **4**. The Plate Handler arm moves over the input stack.



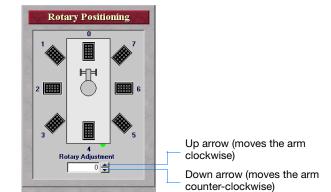
- **3.** Using the Vertical Positioning commands, lower the Plate Handler arm until it is just above the stack. The Vertical Positioning commands provide five ways to move the Plate Handler arm:
  - Move the slider for large increments.
  - Click inside the slider bar to move the arm in 250-step increments.
  - Click the lower arrow on the bar to move the arm in 50-step increments.
  - Click the up or down arrows in the Vertical Adjustment text box to move the arm in 1-step increments.
  - Click the **Vertical Adjustment** text box, enter a value, then press **Enter** to move the arm into a specific location.



- **4.** Using the Rotary Adjustment arrows, adjust the rotational position of the gripper so that it is centered over the input stack and will not contact the sides when lowered.
  - To move the Plate Handler arm clockwise, click the up arrow.
  - To move the Plate Handler arm counterclockwise, click the down arrow.
- 5. Using the Vertical Positioning commands, carefully lower the Plate Handler arm into the stack. Change the Rotary Adjustment value as needed to center the gripper inside the stack.
- 6. After the gripper is centered inside the stack, click to lower the Plate Handler arm on the plate.
- **7.** Confirm the following:
  - The plate is in the middle of the gripper span
  - The plate sensor switch is contacting the plate
  - The gripper and plate do not contact the side of the stack
- **8.** Click Close Gripper . The gripper grips the plate between its fingers.
- 9. Select to Vertical Home. The Plate Handler raises the arm to its highest position. If the plate contacts the sides of the stack, readjust the rotary position of the Plate Handler arm until the plate moves freely in the stack.

**Note:** Contact between the plate and the stack or all stacks may be unavoidable. However, try to minimize the contact as much as possible.

**10.** Using the Vertical Positioning commands, raise and lower the Plate Handler arm several times to check the alignment.





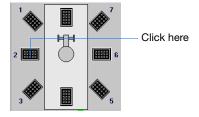
- **11.** Lower the Plate Handler arm to the bottom of the plate stack, click Rotary Offset, then click Yes. The software records the rotary position for the Zymark position 4 (input stack 1).
- **12.** Click Open Gripper. The gripper releases the plate.

# Aligning the Plate Handler to the Instrument Tray

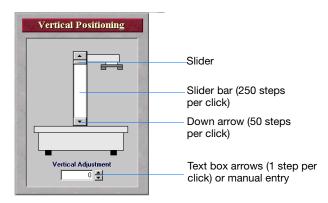
The next step is to align the Plate Handler arm to the instrument tray (Zymark position 2). This alignment will ensure a smooth exchange between the Plate Handler arm and the instrument tray during instrument operation.

- **1.** If not already present, place an empty plate into input stack 1 (Zymark position 4) and pick it up with the Plate Handler arm:
  - a. In the Zymark Twister Software, click position 4.

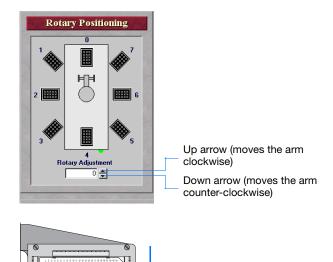
  - c. Click Close Gripper
- **2.** Click position **2**. The Plate Handler arm moves over the instrument tray.



**3.** Using the Vertical Positioning commands, lower the Plate Handler arm until it is approximately 1 cm above the instrument tray.

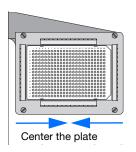


**4.** Using the Rotary Adjustment arrows, center the gripper and plate along the Y-axis of the instrument tray.

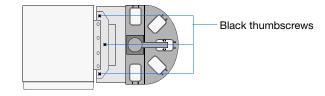


Center the plate

**5.** Center the gripper and plate along the X-axis of the instrument tray by sliding the Plate Handler and base towards or away from the 7900HT Fast System.



- **6.** Again using the Rotary Adjustment arrows, center the gripper and plate along the Y-axis of the instrument tray, as shown in step 4 above.
- 7. Using the Vertical Positioning commands, carefully lower the Plate Handler arm onto the instrument tray and confirm that the plate rests squarely inside it.
- **8.** Tighten the three black thumbscrews on the platform connecting the 7900HT Fast System and the Plate Handler.



- **9.** Release the plate from the Plate Handler arm:
  - a. Click Open Gripper
  - b. Click to Vertical Home .

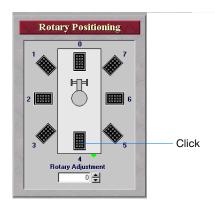


- **10.** Click . The Plate Handler arm lowers onto the plate.
- **11.** Save the rotary and vertical offset information:
  - a. Click Rotary Offset, then click Yes. The software records the rotary position for the plate drawer (Zymark position 2).
  - b. Click vertical Offset, then click vertical position for the plate drawer.

## Rechecking the Input Stack 1

Now that the positions of the Plate Handler and instrument are fixed, the Plate Handler stacks can be aligned and the positional values recorded.

- **1.** If not already present, place an empty plate into input stack 1 (Zymark position 4).
- **2.** In the Zymark Twister Software, click position **4**. The Plate Handler arm moves over the input stack.



- **3.** Using the Vertical Positioning commands, lower the Plate Handler arm until it is 1 cm above the stack and verify that it is centered on the stack. If necessary, center the stack using the Rotary Adjustment arrows.
- **4.** Carefully lower the Plate Handler arm into the stack. Center the gripper as it moves down the stack by adjusting the Rotary Adjustment arrows if needed.
- **5.** After the Plate Handler arm is centered inside the stack, click Find Plate . The Plate Handler arm lowers upon the plate.

- **6.** Confirm the following:
  - The plate is in the middle of the gripper span.
  - The plate sensor switch is contacting the plate.
  - The gripper does not contact the side of the stack.
- 7. Click 

  □ Close Gripper
- **8.** Click to Vertical Home. The Plate Handler raises the arm to its highest position. If the plate contacts the sides of the stack, re-adjust the rotary position of the Plate Handler arm until the plate moves freely inside the stack.

**Note:** Contact between the plate and the stack or all stacks may be unavoidable. However, try to minimize the contact as much as possible.

- 9. Click Rotary Offset, then click Yes. The software re-records the rotary position for input stack 1 (Zymark position 4).
- **10.** While holding the plate, click per Gripper and remove the plate.

# Defining the Bottom of the Stack

The Automation Controller Software requires a bottom position value for all stacks. This value is used to prevent the Plate Handler arm from colliding or grinding as it moves to the bottom of each stack.

- 1. Remove all plates from the instrument and the Plate Handler arm.
- **2.** Place an empty plate into the output stack (Zymark position 0).
- **3.** In the Zymark Twister Software, click position **0**. The Plate Handler arm moves over the output stack.
- **4.** Using the Vertical Positioning commands, lower the Plate Handler arm until it is just above the stack.
- **5.** Using the Rotary Adjustment arrows, adjust the rotational position of the gripper so that it is centered over the input stack and will not contact the sides when lowered.
- **6.** Using the Vertical Positioning commands, carefully lower the Plate Handler arm into the stack. Change the Rotary Adjustment value as needed to center the gripper inside the stack.
- 7. After the gripper is centered inside the stack, click . The Plate Handler arm lowers upon the plate.



- **8.** Confirm the following:
  - The plate is in the middle of the gripper span
  - The plate sensor switch is contacting the plate
  - The gripper does not contact the side of the stack
- 9. Click Gripper |. The gripper grips the plate between its fingers.
- **10.** Select to Vertical Home. The Plate Handler raises the arm to its highest position. If the plate contacts the sides of the stack, readjust the rotary position of the Plate Handler arm until the plate moves freely in the stack.

**Note:** Contact between the plate and the stack may be unavoidable. However, try to minimize the contact as much as possible.

- **11.** Using the Vertical Positioning commands, raise and lower Plate Handler arm several times to check the alignment.
- **12.** Lower the Plate Handler arm, click Rotary Offset, then click Yes. The software records the rotary position for position 0 (the output stack).
- 13. Click the tvertical Home
- **14.** While holding the plate, click open Gripper and remove the plate.
- **15.** In the Vertical Adjustment field, enter **–3200**, then press **Enter**. The Plate Handler lowers the arm to a position near the base of the output stack.
- **16.** Carefully lower the Plate Handler arm until it is approximately 1–2 mm from the bottom of the stack.
- **17.** Click vertical Offset, click vertical position for position 0 (the output stack).
- **18.** Click to Vertical Home. The Plate Handler raises the Plate Handler arm to its highest position.
- **19.** In the Vertical Adjustment field, enter the Vertical Offset value determined in step 17, and press **Enter**. The Plate Handler lowers the Plate Handler arm to a Vertical Offset position.
- **20.** If necessary, readjust the **Vertical Offset** value and repeat steps 18 through 19 until satisfied with the setting.

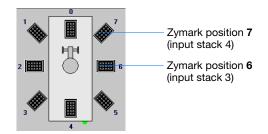
- **1.** Place an empty plate into input stack 2 (Zymark position 5).
- **2.** In the Zymark Twister Software, click position **5**. The Plate Handler arm moves over the input stack.
- **3.** Using the Vertical Positioning commands, lower the Plate Handler arm until it is approximately 1 cm above the stack, then center it using the Rotary Adjustment arrows.
- **4.** Carefully lower the Plate Handler arm into the stack. Center the gripper as it moves down the stack by adjusting the Rotary Adjustment arrows as needed.
- **5.** After the Plate Handler arm is centered inside the stack, click Plate Handler arm lowers upon the plate.
- **6.** Confirm the following:
  - The plate is in the middle of the gripper span.
  - The plate sensor switch is contacting the plate.
  - The gripper does not contact the side of the stack.
- 7. Click Close Gripper
- 8. Click to Vertical Home. The Plate Handler arm raises to its highest position. If the plate contacts the sides of the stack, readjust the rotary position of the Plate Handler arm until the plate moves freely in the stack.

**Note:** Contact between the plate and the stack may be unavoidable. However, try to minimize the contact as much as possible.

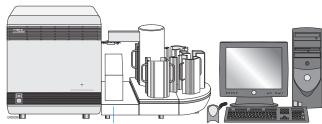
- **9.** Using the **Vertical Positioning** commands, raise and lower Plate Handler arm several times to check the alignment.
- 10. Click Rotary Offset, and click Yes. The software records the rotary position for the Zymark position 5 (input stack 2).



**11.** Repeat steps 1 through 10 for input stacks 3 and 4 to define Rotary Offset values for the remaining positions 6 and 7.



- **12.** Exit the Zymark Twister Software:
  - a. Click Main Menu
  - **b.** Click **Quit Application** to close the software.
- **13.** Replace the covers for the fixed-position bar code reader and the underlying platform (removed in step 1 on page 99).



Fixed-position bar code reader and underlying platform covers

# Confirm the Zymark Twister Software Is Closed

A defect in the Zymark Twister Software can cause portions of the program to persist in memory even after the software has been closed. Because the Zymark Twister Software conflicts with the SDS software, the residual elements of the Zymark Twister Software must be closed inside the Windows operating system Task Manager before continuing.

- **1.** Press the **Crtl** + **Alt** + **Del** keys in unison to open the Windows operating system Security dialog box.
- **2.** Click **Task Manager** to open the Windows operating system Task Manager dialog box.
- **3.** Click the **Applications** tab.

-N	Л	$\overline{}$	+	$\overline{}$	c
Ш	И	U	ι	U	c

- **4.** Confirm that the software has closed by looking for the Zymark Twister Software entry in the **Task** list.
  - If the software is completely closed, there is no entry for the Zymark Twister Software.
  - If the software is still running, click the Zymark Twister Software entry to select it, then click **End Task**.
- **5.** Close the Task Manager dialog box.

## Aligning the Fixed-Position Bar Code Reader

### Two Bar Code Readers

Two bar code readers are available for the 7900HT Fast System:

- CLV Fixed-Position Bar Code Reader
- LD Fixed-Position Bar Code Reader

This section provides procedures for aligning both bar code readers.

#### When to Perform

The fixed-position bar code reader must be set so that it automatically scans the plate's bar code as the plate is placed into the instrument tray by the Plate Handler.

Align the fixed-position bar code reader:

• Whenever the Automation Accessory platform is moved from its calibrated position (for example, during Service visits)

## Before You Begin

## Time Required

20 minutes

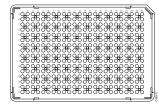
## Materials Required

You need the materials listed below to perform this procedure.

Appropriate plate type for your current sample block; one of the following:

- Standard 384-well reaction plate
- Standard 96-well reaction plate
- TaqMan® Low Density Array

**Note:** The Automation Accessory should not be used for running Fast 96-well reaction plates.



Pictured: Standard 384-well reaction plate

#### Workflow

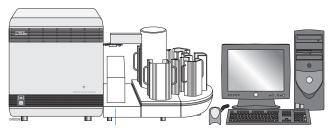
#### To align the fixed-position bar code reader:

- **1.** Obtain the tools and materials required (shown above).
- **2.** Prepare the 7900HT instrument (see page 111).
- **3.** Align the fixed-position bar code reader. Be sure to follow the right alignment procedure for your bar code reader:
  - "Aligning the CLV Fixed-Position Bar Code Reader" on page 112, or
  - "Aligning the LD Fixed-Position Bar Code Reader" on page 114

Notes		

## Preparing the Instrument for the Alignment

**1.** Remove the cover for the fixed-position bar code reader.



Fixed-position bar code reader cover

- **2.** Power on the 7900HT Fast System and the computer.
- **3.** Move the instrument tray to the OUT position:
  - a. Double-click on the desktop (or select Start > All Programs > Applied
     Biosystems > SDS 2.3 > SDS Automation
     Controller 2.3) to start the Automation
     Controller Software.

Note: If an error dialog box appears reading, Machine calibration values are not valid. Please refer to documentation for calibration process, click \_\_\_\_ok\_\_\_.

- **b.** Click **Open/Close Door**. The 7900HT Fast System moves the instrument tray to the OUT position.
- **c.** Select **File** > **Exit** to close the Automation Controller Software.

#### **4.** Continue with:

- "Aligning the CLV Fixed-Position Bar Code Reader" on page 112, *or*
- "Aligning the LD Fixed-Position Bar Code Reader" on page 114

## Aligning the CLV Fixed-Position Bar Code Reader

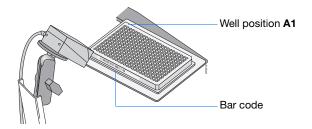
**Note:** This procedure is for the CLV Fixed-Position Bar Code Reader. If you are aligning the LD Fixed-Position Bar Code Reader, see page 114.

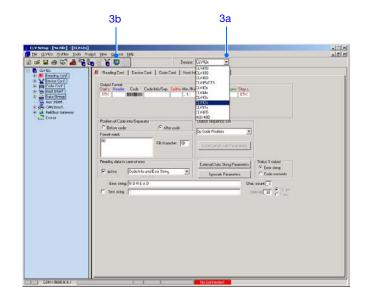
**IMPORTANT!** The instrument tray must be in the OUT position to align the bar code reader.

**1.** Place a plate with bar code onto the instrument tray.

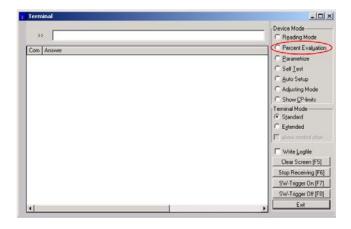
**IMPORTANT!** Orient the plate so that well **A1** aligns to the **A1** position of the instrument tray and that the bar code faces the fixed-position bar code reader.

- 2. Select Start > All Programs > CLV Setup v4.1 to start the CLV software.
- **3.** In the CLV Setup window:
  - a. Click the **Device** drop-down list and select **CLV42x**.
  - **b.** Click **(Terminal)** to open the Terminal dialog box.





4. In the Terminal dialog box, select **Percent** Evaluation.

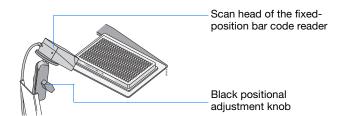


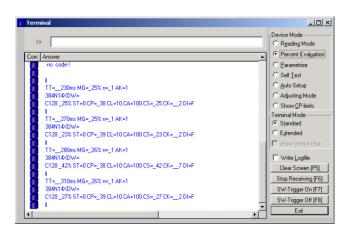
- **5.** Loosen the black positional adjustment knob on the fixed-position bar code reader, and position the scan head of the reader as far as possible from the plate while maintaining the orientation towards the bar code on the plate.
- **6.** While watching the Terminal dialog box, slowly adjust the orientation of the fixed-position bar code reader until the Terminal dialog box reads more than 20%.

**Note:** The Percent Evaluation depends on how many lines hit the bar code label.

**Note:** It may be helpful to briefly place a sheet of white paper in front of the plate bar code to view the area scanned by the laser.

- **7.** When satisfied with the alignment, tighten the black positional adjustment knob on the fixed-position bar code reader.
- **8.** Click **Exit** to close the Terminal dialog box, then close the CLV software.
- **9.** Replace the cover for the fixed-position bar code reader (removed in step 1 on page 111).





## Aligning the LD Fixed-Position Bar Code Reader

**Note:** This procedure is for the LD Fixed-Position Bar Code Reader. If you are aligning the CLV Fixed-Position Bar Code Reader, see page 112.

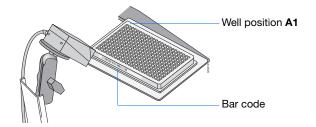
**IMPORTANT!** The instrument tray must be in the OUT position to align the bar code reader.

**1.** Place a plate with bar code onto the instrument tray.

**IMPORTANT!** Orient the plate so that well **A1** aligns to the **A1** position of the instrument tray and that the bar code faces the fixed-position bar code reader.

- 2. Select Start > All Programs > PSC Laser Data > LDHOST to start the LDHOST software.
- **3.** Establish communication with the fixed-position bar code reader:
  - a. Click [2] (**Edit**) to open the Edit Configuration dialog box.
  - b. Click (Terminal) to open the Terminal dialog box.
  - c. In the Device Control dialog box, click (Connect to Device). The Terminal dialog box displays the fixed-position bar code reader response.
  - d. Click or to close the information message.

The LD Host program communicates with the bar code reader and updates the Edit Configuration dialog box with the current configuration settings.



## **4.** Configure the software for the alignment:

 a. In the bottom of the Edit Configuration dialog box, locate and select the Op. Modes tab.

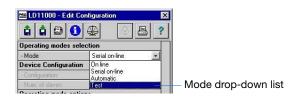
**Note:** You may need to use the arrows located in the bottom of the dialog box to locate the Op. Modes tab.

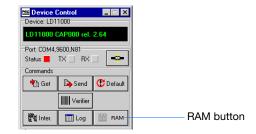
- b. In the Operating modes selection pane of the Edit Configuration dialog box, click the Mode drop-down list and select Test.
- c. In the Device Control dialog box, click RAM to toggle to EEPROM mode, then click Send.
- **d.** In the Confirm dialog box, click **YES** to save to EEPROM.

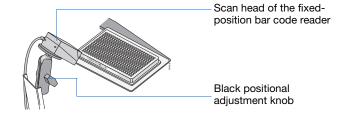
The fixed-position bar code reader begins a continuous repeating scan of the bar code. The software updates the Terminal dialog box every 0.5 sec indicating the percentage of accurate reads completed during the 0.5 sec interval.

5. Loosen the black positional adjustment knob on the fixed-position bar code reader, and position the scan head of the reader as far as possible from the plate while maintaining the orientation towards the bar code on the plate.











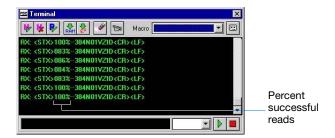
**6.** While watching the Terminal dialog box, slowly adjust the orientation of the fixed-position bar code reader until the percent successful reading displays the highest number possible.

**Note:** It may be helpful to briefly place a sheet of white paper in front of the plate bar code to view the area scanned by the laser.

- **7.** When satisfied with the alignment, tighten the black positional adjustment knob on the fixed-position bar code reader.
- **8.** Restore the fixed-position bar code reader to normal operation:
  - a. In the Operating modes selection pane of the Edit Configuration dialog box, click the Mode drop-down list and select Serial on Line.
  - b. In the Device Control dialog box, confirm that **EEPROM** is still selected, then click **Send**
  - **c.** In the New Decision dialog box, click **YES** to save to EEPROM.

The bar code reader stops scanning the plate bar code and resumes normal operation.

- **9.** Click **(Exit)** to close the LDHOST software.
- **10.** Replace the cover for the fixed-position bar code reader (removed in step 1 on page 111).





Mode drop-down list

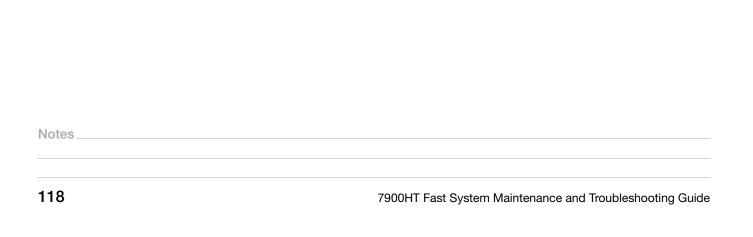
## **Troubleshooting the Automation Accessory**

# Troubleshooting Table

Table 4-2 Troubleshooting the Automation Accessory

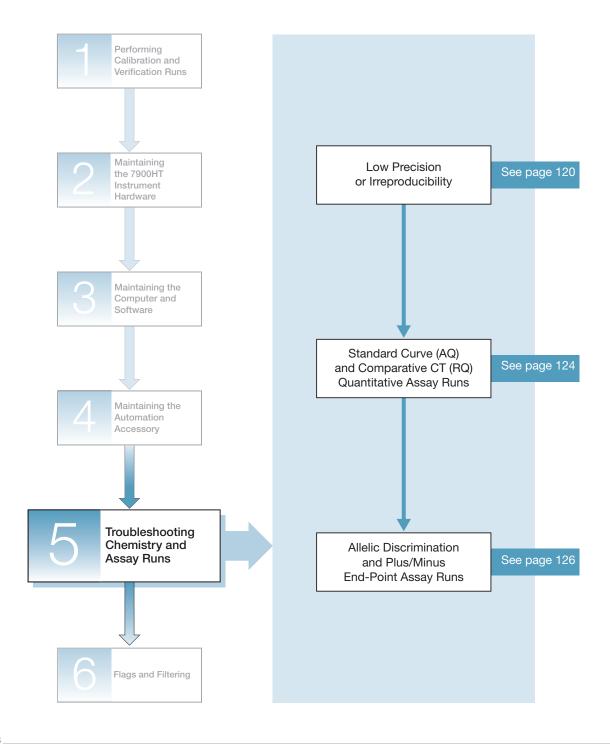
Observation	Possible Cause	Recommended Action		
Plate Handler emits	Vertical offset too low	Re-align the Plate Handler as		
grinding noise when picking up or putting down plates	Plate detector switch set too high	explained in "Aligning the Plate Handler" on page 97.		
Plate Handler arm contacts racks when retrieving or stacking plates	Plate Handler rotary offset is incorrect or vertical offset is too low			
The Plate Handler arm releases plates awkwardly into the plate racks				
Reaction plates tip or tilt when placed into the instrument tray by the Plate Handler arm				
Plate Handler fails to sense or grasp plates	Plate sensor switch not adjusted properly	Adjust the plate sensor switch as explained in "Adjusting the Plate Sensor Switch" on page 92.		
	Gripper pads on the fingers of the Plate Handler arm are worn or dirty	Change the gripper pads as explained in "Cleaning and Replacing Gripper Finger Pads" on page 90.		
Plates stick to the gripper fingers of the Plate Handler arm	Gripper pads are worn or dirty	Change the gripper pads as explained in "Cleaning and Replacing Gripper Finger Pads" on page 90.		
Plate Handler does not restack plates in original locations	Restack when finished option not selected	Configure the Automation Controller Software to restack the plates. See the Sequence Detection Systems Software version 2.3 Online Help (SDS Online Help).		
Fixed-position bar code reader not reading plate	Bar code reader is misaligned	Re-align the fixed-position bar code reader as explained in		
bar codes	Bar code reader is broken	"Aligning the Fixed-Position Bar Code Reader" on page 110.		

# Chapter 4 Maintaining the Automation Accessory Troubleshooting the Automation Accessory



# 5

# Troubleshooting Chemistry and Assay Runs



## Low Precision or Irreproducibility

#### Overview

There are various reasons why an assay run with the Applied Biosystems 7900HT Fast Real-Time PCR System can have less than optimal precision. Factors that can affect precision are described in detail below.

Improper Threshold Setting	120
Imprecise Pipetting	121
Non-Optimized Chemistry	121
Incomplete Mixing	121
Air Bubbles	121
Splashing PCR Reagents	121
Drops	122
Writing on the Reaction Plates	122
Fluorescent Contamination on the Plates	122
Errors	122
Contaminated Sample Block	122
Improper or Damaged Plastics	123
Low Copy Templates	123
Use of Non-Applied Biosystems PCR Reagents	123

## Improper Threshold Setting

The key to high-precision quantitative PCR is accurate detection of the exponential (geometric) phase. The Applied Biosystems 7900HT Fast Real-Time PCR System typically delivers sufficient sensitivity so that at least three cycles of the geometric phase are visible, assuming reasonably optimized PCR conditions. The SDS software calculates a fixed signal intensity, called a threshold, that each signal generated from PCR amplification must reach before it is recognized by the software as actual amplification. The calculated threshold is an approximation, and should be examined and modified as needed.

#### Modifying the Threshold

In a real-time document of the SDS software, the threshold can be modified via the Amplification Plot view following analysis of the run data. See the *Sequence Detection Systems Software version 2.3 Online Help (SDS Online Help)* for more information.

Notes		

# Imprecise Pipetting

The calculated quantities of target nucleic acid are directly affected by how precisely the template or reagent volumes are added to the reaction mixes.

#### **Using Master Reaction Mixes**

Applied Biosystems highly recommends using a master reaction mixture. Mix all common components (including the same template) to a set of reactions together, then dispense the mix to the replicate wells of the plate. When making each master reaction mix, add 5–10% additional volume to compensate for pipetting losses.

#### **Using Pipettors**

Pipetting precision is also improved by:

- · Calibrating and servicing the pipettors regularly
- Pipetting larger volumes
- Reducing the number of pipetting steps whenever possible
- Increasing the consistency of the pipetting method

Consult the manufacturer about the correct method of dispensing liquid volumes accurately from the pipettor. For example, some pipettors are designed to deliver the designated volume at the first plunger stop, so 'blowing out' the residue may cause error. Also, before using a new pipettor tip to serially dispense a master reaction mix, wet the tip once by drawing up some of the master reaction mix and dispensing it back into the mix again.

# Non-Optimized Chemistry

Chemistries that have not been optimized may be susceptible to inconsistencies. To maximize precision and reaction efficiency, optimize the primer and probe concentrations of each individual assay used. Refer to the *TaqMan® Universal PCR Master Mix Protocol* (PN 4304449) for specific information about optimizing probe and primer concentrations for TaqMan-related chemistries.

**Note:** TaqMan<sup>®</sup> Gene Expression Assays, TaqMan<sup>®</sup> SNP Genotyping Assays, and Custom TaqMan<sup>®</sup> SNP Genotyping Assays do not require optimization.

# Incomplete Mixing

For maximum precision, the PCR master reaction mix must be uniform. After adding all reaction components to the mix, vortex the mix for 4–5 seconds before aliquoting it to the wells of the plate. Also vortex any dilutions performed during the assay.

#### Air Bubbles

Air bubbles in the wells can refract and distort the fluorescent signals. Ideally, the reagents would be applied to the wells using a pipetting technique that does not form air bubbles. However, if a plate does contain air bubbles, they can usually be removed by swinging, tapping, or briefly centrifuging the reaction plate.

# Splashing PCR Reagents

If PCR reagents splash the undersides of the optical adhesive covers, the heat from the lid may bake the liquid to the cover and may distort the signal. If splashing occurs, briefly centrifuge the reaction plate to remove all traces of liquid from the caps.



#### **Drops**

Drops of reagents that cling to the sides of the wells may not contact the thermal cycler sample block and consequently may not amplify. If the drop slides into the mix during PCR, then the amplified products will become diluted and the result will be less than replicate wells that did not have drops. Therefore, carefully monitor the reaction plate as it is being transferred into the thermal cycler or 7900HT Fast System. If you observe any drops, take steps to remove them, such as centrifugation.

## Writing on the Reaction Plates

Do not write on any surface of the optical plates or the optical adhesive covers. The fluorescent properties of the ink can potentially affect the fluorescence emission from the plate and alter the results. Instead, note the contents of each well on a sheet of paper, or on a printout of the sample setup.

## Fluorescent Contamination on the Plates

Many compounds found in laboratories are fluorescent. If they come into contact with certain optical surfaces, such as the optical adhesive covers, the fluorescent results may be affected. For example, it has been noted that the powder used to lubricate the insides of plastic gloves often contains fluorescent compounds. Use only powder-free gloves and do not needlessly touch the reaction plates or optical adhesive seals.

#### **Errors**

Human errors from time to time are inevitable, such as pipetting into the wrong well, or making a dilution mistake.

Human error can be reduced in the following ways:

- Perform the assay in a systematic fashion. For example, the pattern of sample positions should be simple (avoid putting gaps in the rows).
- When pipetting the master reaction mix, look directly down into the reaction plate so that you can verify the transfer of the solution.
- If adding a small-volume reagent, such as template, place the drop of liquid on the side of the well. Briefly tap or centrifuge the plate afterwards to bring the droplet down into the well.
- After all pipetting is complete, visually inspect all the wells to confirm the presence of the reagent drops. Tapping or centrifuging the reaction plate will cause all the drops to slide down into the wells simultaneously.
- When making serial dilutions, be sure to change the pipet tip after each dilution step.
- Visually inspect the liquid volumes being pipetted to verify that the volume is approximately correct. A common mistake is using the wrong pipettor volume setting (such as setting 20 µL instead of 2.0 µL).
- Visually inspect the volumes of the completed reactions, looking for any wells that have volumes that do not match those of the other wells.

## Contaminated Sample Block

Any material contaminating the sample block can affect the results. For example, mineral oil reduces thermal transfer. Residue from writing on reaction plates darkens the wells, absorbing light.

Notes		

The sample blocks should be periodically inspected for cleanliness. Sample block contamination can be visualized by running a background plate and inspecting the resulting background signal for aberrant peaks above 4000 FSU (see page 15). See page 65 for instructions on decontaminating the sample block.

## Improper or Damaged Plastics

Applied Biosystems recommends that you use Applied Biosystems optical plates, optical adhesive covers, and optical flat caps with the 7900HT Fast System. The plastics that comprise the optical parts undergo special testing for the absence of fluorescent impurities. Optical plates are frosted to improve the degree and precision of light reflection. Bent, creased, or damaged plastics may adversely affect the transmission of fluorescent signal or prevent proper sealing of a well resulting in evaporation, change in sample volume, and altered PCR chemistry. Make sure to use the correct plastics and visually inspect each reaction plate before use.

**Note:** See Appendix C, "Parts and Consumables," on page 157 for a list of compatible consumables and reagents.

## Low Copy Templates

When amplifying samples that contain very low quantities of nucleic acid (generally less than 100 molecules), expect lowered precision due to the Poisson distribution and biochemical effects related to binding probabilities. Low copy templates are also more susceptible to losses due to non-specific adhesion to plastic wells, pipettor tips, etc. The addition of carrier to the sample, such as yeast tRNA or glycogen, can help prevent these losses, increasing the precision and sensitivity of the assay.

## Use of Non-Applied Biosystems PCR Reagents

The Applied Biosystems buffer contains an internal passive reference molecule  $(ROX^{TM} dye)$ , which acts as a normalization factor for fluorescent emissions detected in the samples.

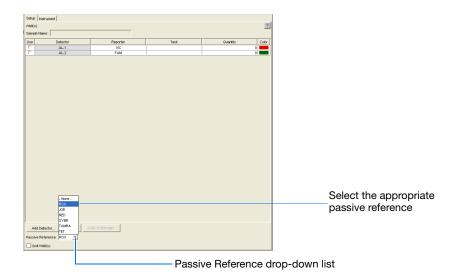
**IMPORTANT!** Non-Applied Biosystems PCR buffers may not contain the ROX passive reference dye. If running non-Applied Biosystems chemistry, be sure to set the passive reference for your experiment as explained below.

#### Setting the Passive Reference

- 1. Open the appropriate plate document in the SDS software.
- 2. Select the **Setup** tab.
- 3. From the Passive Reference drop-down list, select the appropriate passive reference dye.

- N	ь	$\overline{}$	4.	~
1	A (	U	U	E:





# Standard Curve (AQ) and Comparative $C_T$ (RQ) Quantitative Assay Runs

## Troubleshooting Analyzed Data

When faced with irregular data, you can use the SDS software to diagnose some chemistry- and instrument-related problems. The following table contains a summary of checks for verifying the integrity of your run data and to help you begin troubleshooting potential problems.

#### Raw Data Plot

The Raw Data Plot displays the raw reporter fluorescence signal (not normalized) for the selected wells during each cycle of the real-time PCR.

#### What to look for:

- **Signal tightness and uniformity** Do the raw spectra signals from replicate groups and controls exhibit similar spectral 'profiles'? If not, the plate or sample block could be contaminated.
- Characteristic signal shape Do the samples peak at the expected wavelengths? For example, samples containing only FAM<sup>™</sup> dye-labeled TaqMan<sup>®</sup> probes should not produce raw fluorescence in the wavelength of a VIC<sup>®</sup> dye component. A signal present in wells that do not contain the dye could indicate that the sample, master mix, or well contains contaminants.
- Characteristic signal growth As you drag the bar through the PCR cycles, do you observe growth as expected? Absent growth curves may indicate a pipetting error (for example, well lacks template) or no amplification.
- **Signal Plateaus** Do any of the signals plateau? Signal plateaus or saturation can be an indication that a well contains too much template or fluorescent signal.

Notes			

## Multicomponent Plot

The Multicomponent Plot displays a plot of normalized multicomponent data from a single well of a real-time run. The plot displays the component dye signals that contribute to the composite signal for the well.

#### What to look for:

- Correct dyes displayed Does the plot display all dyes as expected? The presence of an unexpected dye may be the result of an error in detector setup, such as assigning the wrong reporter or quencher dye.
- ROX dye fluorescence level Does the ROX dye signal fluoresce below the reporter dyes? If not, the lack of reporter fluorescence may be caused by an absence of probe in the well (a pipetting error).
- Background fluorescence Do all dyes fluoresce above the background? The Background signal is a measure of ambient fluorescence. If a dye fails to fluoresce above the background, it is a strong indication that the well is missing probes labeled with the dye (well does not contain probe, PCR master mix, or both).
- MSE Level The MSE (mean squared error) is a mathematical representation of how accurately the multicomponented data fits the raw data. The higher the MSE value, the greater the deviation between the multicomponented data and the raw data.

## **Amplification Plot**

The Amplification plot displays data from real-time runs after signal normalization and Multicomponent analysis. It contains the tools for setting the baseline and threshold cycle  $(C_T)$  values for the run.

#### What to look for:

• Correct baseline and threshold settings – Are the baseline and threshold values set correctly?

Identify the components of the amplification curve and set the baseline so that the amplification curve growth begins at a cycle number that is greater than the highest baseline number.

Identify the components of the amplification curve and set the threshold so that it is:

- Above the background
- Below the plateaued and linear regions
- Within in the geometric phase of the amplification curve

**IMPORTANT!** After analysis, you must verify that the baseline and threshold were called correctly for each well by clicking on the well in the Plate Grid and viewing the resulting plots. For more information about manually adjusting the baseline and threshold settings, refer to the *SDS Online Help*.

• **Irregular amplification** – Do all samples appear to have amplified normally? The three phases of the amplification curve should be clearly visible in each signal.

Notes			

Outlying amplification – When the run data is viewed in the C<sub>T</sub> vs. Well Position
plot, do replicate wells amplify comparably? Wells producing C<sub>T</sub> values that differ
significantly from the average for the associated replicate wells may be considered
outliers

If a plate produces non-uniformity between replicates, some samples on the plate could have evaporated. Check the seal of the optical adhesive cover for leaks.

## Allelic Discrimination and Plus/Minus End-Point Assay Runs

## Troubleshooting Analyzed Data

When faced with irregular data, you can use the SDS software to diagnose some chemistry- and instrument-related problems. The following table contains a summary of checks for verifying the integrity of your run data and to help you begin troubleshooting potential problems.

#### Raw Data

The Raw Data Plot displays the raw reporter fluorescence signal (not normalized) for the selected wells during each cycle of the PCR.

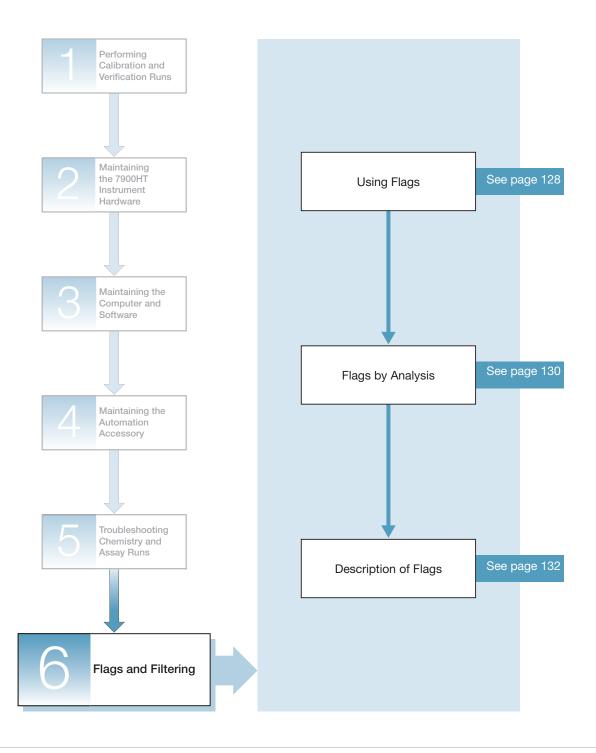
#### What to look for...

- **Signal tightness and uniformity** Do the raw spectra signals from replicate groups and controls exhibit similar spectral 'profiles'? If not, the plate or sample block could be contaminated.
- Characteristic signal shape Do the samples peak at the expected wavelengths? For example, samples containing only FAM dye-labeled TaqMan probes should not produce raw fluorescence in the peak wavelength of the VIC dye component. A signal present in wells that do not contain the dye could indicate that the sample, master mix, or well contains contaminants.
- **Signal Plateaus** Do any of the signals plateau? Signal plateaus or saturation can be an indication that a well contains too much template or fluorescent signal.

Votes			



# Flags and Filtering



## **Using Flags**

## Flags and Troubleshooting

Version 2.3 of the Sequence Detection Systems (SDS) Software features a system of flags that provide the basis for algorithmically identifying and eliminating problematic data from plate document and RQ study analysis. The flags are application-specific metrics, where each evaluates the data for a specific quality that is consistent with a problem common to a supported analysis. The system is designed to allow you to automatically exclude questionable data, alert you to potential problems, and provide you with a starting point for investigation.

## **About Flags**

Each flag is the result of a unique algorithmic test that evaluates a specific property of the plate document data. The software performs the tests in a specific sequence and completes the analysis of the plate document even if one or more wells is flagged. Every flag yields a pass/fail result that indicates the success of the associated test. If a flag is enabled, the software indicates the status of the test in the appropriate column of the Results Table, as follows:

- A **v** for a well that is flagged (Fail)
- Blank for a well that is not flagged (Pass)

In the SDS Software, you can automatically flag results data to meet specified criteria. Flags are assay specific and are user configured or automatically assigned.

- Automatically assigned flags are not user configurable. These flags are displayed in the Results table and the QC Summary tab.
- User configured flags are set/defined by the user. These flags are also displayed in the Results table and QC Summary tab.

#### Flag Symbols

Following an analysis, the software summarizes the results of the flag metrics for each sample by displaying one of four symbols. These symbols appear in the Flag column of the Results Table and in the Plate Grid.

Symbol	Definition
	The sample passed all applicable flag tests (Pass).
2	One or more possible problems exists for the associated sample (Fail).  Note: The symbol displays the number of flags the sample failed.
1	Omitted Well: Well omitted automatically by the algorithm  Note: The symbol displays the number of flags the sample failed.
$\boxtimes$	Omitted Well: Well omitted manually by user

Note:	Applied	Biosystems	recommends	examining	all sampl	es that are	flagged.
-------	---------	------------	------------	-----------	-----------	-------------	----------

Notes\_\_\_\_\_

## **Customizing Flags**

Most flags are a measurement of a specific aspect of the sample data. A flag is produced when the measurement exceeds a threshold, some of which you can customize to refine your analyses. In these cases, the user-defined parameters for the flags appear in the Analysis Settings dialog box for the applicable analysis type. See "Configuring Flags for Use" on page 129 for more information on entering custom thresholds.

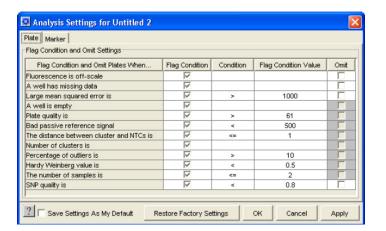
## Notes for Using Flags

- Flags appear in the results tab, the plate grid, and the QC Summary tab.
- Flags are warnings. The software does not stop the analysis if one or more samples fail the flag tests.
- Applied Biosystems recommends examining flagged results ( ).
- Holding the cursor over a column header in the Results Table displays a tooltip that lists the full name of the column (the default names are often acronyms).

# Configuring Flags for Use

While viewing the desired plate document or relative quantitation (RQ) study:

- 1. Click to display the Analysis Settings dialog box.
- **2.** In the Analysis Settings dialog box, select the **Plate** tab (or select the **Study** tab if you are configuring the settings for an RQ study).
- **3.** Using the table on page 130 as a reference, do the following for each row in the Flag Condition and Omit Settings table:
  - **a.** If you want the software to display the results of the flag in the Results Table, select the check box in the Flag Condition column.
  - **b.** Do not change the qualifying statement in the Condition column.
  - **c.** In the Flag Condition Value column, enter a value for the software to apply as the threshold or limit for the associated flag.
  - **d.** If you want the software to omit the wells that fail the flag from the analysis, select the check box in the Omit column.



Notes

6

- **4.** If you want the software to apply the current settings to all future plate documents of the same analysis type, select **Save Settings As My Default**.
- **5.** Click **Apply** to reanalyze the plate document or study without closing the Analysis Settings dialog box, or click **OK**.

## Flags By Analysis

**Note:** The plus/minus plate documents do not feature a system of flags.

## Flags for SDS Software Version 2.3

Flags		rable lute ation ic		ic nation	ic nation inus	Relative Quantitation		See
Abb.	Name	User Configurable	User Configurable Absolute Quantitation	Absolute Quantitation Allelic Discrimination	Plus/Minus	Plate Docu- ment	Study	Page
BAF	Baselining algorithm failed		~				<b>'</b>	132
BPR	Bad passive reference signal	~	~	~		<b>V</b>	<b>'</b>	132
CAF	C <sub>T</sub> calculation algorithm failed		~				<b>'</b>	133
DCN	The distance between cluster and NTCs is	~		~				133
EAF	Exponential region algorithm failed		~				<b>'</b>	133
EW	A well is empty	~	~	~		<b>V</b>	<b>'</b>	134
FOS	Fluorescence is off-scale	~	<b>'</b>	<b>'</b>		~	<b>'</b>	134
GBO	Well is an outlier	~					<b>~</b>	139
HMD	A well has missing data	~	<b>'</b>	<b>'</b>		~	<b>'</b>	136
HNS	A well has a noise spike	<b>'</b>	<b>'</b>			~	<b>'</b>	136
HRN	A well has high relative noise	<b>'</b>	<b>'</b>			~	<b>'</b>	136
HSD	High standard deviation in replicate group	~					~	137
HW	Hardy Weinberg value is	~		~				135
LME	Large mean squared error is	~	~	~		<b>V</b>	~	137
LPL	Laser power is low during the run	<b>'</b>	~			<b>'</b>	<b>✓</b> ‡	138
NAP	Percentage of plate wells not amplified	<b>'</b>	<b>v</b>			~	<b>✓</b> ‡	138

Notes		

## Flags for SDS Software Version 2.3

Flags		rrable ute ation	ute ation	ute ation ic nation	inus	Relative Quantitation		See
Abb.	Name	User Configurable	Absolute Quantitation	Allelic Discrimination	Plus/Minus	Plate Docu- ment	Study	Page
NAW	A well is not amplified	V	~			<b>/</b>	~	138
NOC	Number of clusters is	V		V				139
POU	Percentage of outliers is	V		V				139
SNS	The number of samples is	V		V				139
TAF	Thresholding region algorithm failed		<b>V</b>				<b>'</b>	140

<sup>‡</sup> Visible only when viewing study data using the Plate Centric Table Orientation.

Notes\_\_\_\_

## **Description of Flags**

## **Baselining Algorithm Failed (BAF)**

## Description/ Function

This flag indicates that the algorithm was unable to determine the best fit baseline for the data set.

**Note:** The Exponential Region, Baselining, Thesholding, and  $C_T$  Calculation Algorithms are performed in sequence. If one flag fails, then the SDS software does not compute the flags for the remaining algorithms in the sequence.

## **Troubleshooting**

- **1.** Select the affected sample in the Results Table.
- 2. Select the **Results** tab
- **3.** In the Amplification Plot, review the amplification curves. If necessary, set the baseline and threshold manually for the affected data.

## **Bad Passive Reference (BPR)**

## Description/ Function

This flag indicates that the passive reference signal for the associated sample may:

- · Be below normal levels
- Contain irregular fluctuations (non-uniform signal)

### **Troubleshooting**

- **1.** Confirm the results of the flag test:
  - a. Select the affected sample in the Results Table.
  - b. Select the **Results** tab.
  - **c.** Click to display the Multicomponent Plot.
  - **d.** In the Multicomponent Plot, review the passive reference signal for irregularities.
- **2.** If the passive reference signal does contain inconsistencies, determine the location of the failed well in the Results Table.
- **3.** If the plate or TLDA used to run the sample is available, verify that the well contains fluid.

Fluctuations in the passive reference signal can be caused by:

- An empty or leaking well
- Formation of bubbles in the well

Notes		

## C<sub>T</sub> Calculation Algorithm Failed (CAF)

## Description/ Function

This flag indicates that the software could not calculate a C<sub>T</sub> because of a computational error.

**Note:** The Exponential Region, Baselining, Thesholding, and  $C_T$  Calculation Algorithms are performed in sequence. If one flag fails, then the SDS software does not compute the flags for the remaining algorithms in the sequence.

## **Troubleshooting**

- **1.** Select the affected sample in the Results Table.
- **2.** Select the **Results** tab.
- **3.** In the Amplification Plot, review the amplification curves. If necessary, set the baseline and threshold manually for the affected data.

## Distance Between Cluster and NTCs (DCN)

## Description/ Function

This flag indicates that the associated sample is located close to the no template control (NTC) cluster, and may not have amplified well. If active, the SDS software omits all samples that do not meet condition criteria.

## Troubleshooting

Insufficient signal may be the result of poor PCR or an insufficient number of thermocycles.

## **Exponential Region Algorithm Failed (EAF)**

## Description/ Function

This flag indicates that the software could not identify the exponential region of the amplification curve.

**Note:** The Exponential Region, Baselining, Thesholding, and C<sub>T</sub> Calculation Algorithms are performed in sequence. If one flag fails, then the SDS software does not compute the flags for the remaining algorithms in the sequence.

## **Troubleshooting**

- **1.** Select the affected sample in the Results Table.
- 2. Select the **Results** tab.

**3.** In the Amplification Plot, review the amplification curves. If necessary, set the baseline and threshold manually for the affected data.

## **Empty Well (EW)**

## Description/ Function

This flag indicates that the software has determined the associated well to be empty because the well:

- Does not contain sample or reagents (primers and TaqMan<sup>®</sup> probes).
- Produced very low fluorescence (or none at all)

## **Troubleshooting**

- **1.** Confirm the results of the flag test:
  - **a.** Select the affected sample in the Results Table.
  - b. Select the **Results** tab.
  - **c.** Click to display the Multicomponent Plot.
  - **d.** In the Multicomponent Plot, verify that the signals are below normal level for the plate/study.
- **2.** In the Results Table, determine the location of the failed well.
- **3.** If the plate or TLDA used to run the sample is available, verify that the well contains fluid.

## Fluorescence Off-Scale (FOS)

## Description/ Function

This flag indicates that the associated sample produced fluorescence in one or more cycles that exceeds the maximum detectable range for the 7900HT instrument (CCD saturation).

**Note:** Offscale data can corrupt the results of a run if the offscale data is located at a cycle that interferes with the  $C_T$  calculation.

## **Troubleshooting**

- **1.** Select the affected sample in the Results Table.
- 2. Select the **Results** tab.
- **3.** Click to display the Multicomponent Plot.
- **4.** In the Raw Data and Multicomponent Plots, evaluate the effect of the offscale data. Does the offscale data interfere with the C<sub>T</sub> calculation? If so, consider omitting the well from the analysis.

Notes			

**5.** If possible, determine the source of the offscale result.

A variety of factors can produce offscale data including:

- Sample Overload High concentrations of sample can produce offscale data.
- Contamination A fluorescent contaminant on the exterior a well or within the reaction can produce consistently high signals throughout a run.

## **Outlier Well (GBO)**

## Description/ Function

This flag indicates that the  $C_T$  of the sample deviates significantly from those of the associated replicate group and has been identified as an outlier based on the Grubbs test (also known as the extreme studentized deviate [ESD] method for outlier removal).

**Note:** The algorithm does not remove samples that achieve  $C_T$ s within 0.25 of the mean for the replicate group.

**Note:** If two sample produce identical  $C_T$ s but are identified as outliers, then the software flags both wells.

## **Troubleshooting**

- **1.** In the SDS RQ Manager Software, note the well positions of the outlier and associated replicates.
- 2. Select Table Orientation ➤ Detector Centric.
- **3.** In the upper-left pane, select the detector that the outlier evaluates to view the data for the replicate group.
- **4.** Select the **Amplification Plot** tab, then select **Data**  $\triangleright$  **Ct vs. Well Position** to view the clustering of the  $C_T$ s for the replicate group.
- **5.** Evaluate the outlier. Omit or restore the well as necessary.

## Hardy Weinberg Value (HW)

## Description/ Function

This flag identifies samples that belong to a population that does not match Hardy Weinberg equilibrium.

## Troubleshooting

The samples may trigger this flag as a result of genuine allelic disequilibrium in the associated population or because of several experimental and technical issues.

Notes		

6

## Has Missing Data (HMD)

## Description/ Function

This flag indicates that the associated sample contains one or more cycles for which no spectral data is present.

**Note:** Although rare, samples can contain missing data as the result of an interruption to the 7900HT instrument hardware or software during the run.

## **Troubleshooting**

- **1.** Select the affected sample in the Results Table.
- **2.** Select the **Results** tab.
- **3.** Click me to display the Multicomponent Plot.
- **4.** In the Raw Data and Multicomponent Plots, observe the missing data.

  Does the missing data interfere with the C<sub>T</sub> calculation? If so, consider omitting the well from the analysis.

## Has Noise Spike (HNS)

## Description/ Function

This flag indicates that the amplification curve contains one or more abnormal data points which are inconsistent with the rest of the curve.

### **Troubleshooting**

- **1.** Select the affected sample in the Results Table.
- **2.** Select the **Results** tab.
- **3.** In the Amplification Plot, review the amplification curves for the noise spike. If the noise spike has biased the baselining or C<sub>T</sub> calculations, consider omitting the well from the analysis or set the baseline and threshold manually for the affected data.
- **4.** If no noise spikes are present, change the analysis settings for HNS to be less sensitive.

## **High Relative Noise (HRN)**

## Description/ Function

This flag indicates that the associated sample failed the high relative noise test.

The algorithm determines the magnitude of the noise for each well based on the reporter dye signal. It also estimates the mean and standard deviation of the noise for the entire plate. The flag is evaluated by comparing the noise of the well to the plate, then determining whether or not it exceeds the threshold in the analysis settings.

Notes			

## **Troubleshooting**

- 1. Select the affected sample in the Results Table.
- **2.** Select the **Results** tab.
- **3.** Click me to display the Multicomponent Plot.
- **4.** In the Multicomponent Plot, observe the signal corresponding to the background fluorescence. If necessary, omit the well.
- **5.** Change the analysis settings to increase or decrease sensitivity, as desired.

# High Standard Deviation in Replicate Group (HSD)

## Description/ Function

This flag indicates that the standard deviation of the  $C_T$ s for the replicate group exceeded the criteria defined in the "High standard deviation in a replicate group" row of the Study tab in the Analysis Settings dialog box.

## **Troubleshooting**

- **1.** In the SDS RQ Manager Software, note the well positions within the outlier and associated replicates.
- 2. Select Table Orientation > Detector Centric.
- **3.** In the upper-left pane, select the appropriate detector to view the relevant data for the replicate group.
- **4.** Select the **Amplification Plot** tab, then select **Data**  $\triangleright$  **Ct vs. Well Position** to view the clustering of the  $C_T$ s for the replicate group.
- **5.** Evaluate the  $C_T$  of the flagged well in relation to the replicates. Omit or restore the well as necessary.

## Large Mean Squared Error (LME)

## Description/ Function

This flag indicates that the associated sample failed the Mean Squared Error test. To perform the test, the software compares the maximum value for the mean squared error for the well to the criteria defined in the "Large Mean squared error is" row the Analysis Settings dialog box. If the statement evaluates *True*, then the software flags the well.

## **Troubleshooting**

- **1.** Select the affected sample in the Results Table.
- 2. Select the **Results** tab.
- 3. Click to display the Multicomponent Plot.

Mata	

**4.** In the Multicomponent Plot, observe the MSE curve in the plot.

Factors that can produce a Large Mean Squared Error include:

- Improper Pure Dye Calibration
- Incorrect dyes specified in setup
- · Offscale fluorescence

## Laser Power Low (LPL)

## Description/ Function

This flag indicates that the power supplied to the laser of the 7900HT instrument decreased below the acceptable level at least once during the associated run.

**Note:** Applicable only to runs performed with version 2.3 and later of the SDS Software.

## **Troubleshooting**

Open the log file for the appropriate 7900HT Fast System and confirm the loss of power during the associated run.

**IMPORTANT!** If the loss of laser power is consistent and is observed over multiple runs, an instrument component may be nearing the end of its life and require service or replacement.

# Non-Amplified Plate (NAP)

## Description/ Function

This flag indicates that the percentage of wells on the plate that did not amplify exceeded the threshold defined in the "Percentage of plate wells not amplified" row of the Analysis Settings dialog box.

**Note:** Empty wells are not included in the test.

## Troubleshooting

Review the data of the wells that failed to amplify (flagged as NAW).

# Not Amplified Well (NAW)

## Description/ Function

This flag indicates that the associated sample failed to amplify. To determine this, the algorithm calculates the difference between the signal at the beginning of the run and the signal at the end of the run.

Applied Biosystems recommends using the default setting for the NAW calculation. To customize the setting, perform successive analyses of your data set using different settings and adjust the setting according to the desired sensitivity.

Notes			

## **Troubleshooting**

- 1. Select the affected sample in the Results Table.
- **2.** Select the **Results** tab.
- **3.** Observe the signals in the Amplification Plot. If necessary, omit the well.

## **Number of Clusters (NOC)**

## Description/ Function

This flag indicates that the number of clusters in the data set (excluding the NTC cluster) is outside the threshold range defined in the "Number of clusters is" row of the Analysis Settings dialog box.

## **Troubleshooting**

- **1.** Select the affected sample in the Results Table.
- **2.** Select the **Results** tab.
- **3.** In the Allele Plot, observe the data point in relation to the related cluster. If necessary, omit the well.

## Percentage of Outliers (POU)

## Description/ Function

This flag indicates that the percentage of outliers on the plate exceeds the threshold defined in the "Percentage of outliers is" row of the Analysis Settings dialog box.

# Small Number of Samples in Cluster (SNS)

## Description/ Function

This flag indicates that the number of data points in the cluster is less than the threshold defined in the "The number of samples is" row of the Analysis Settings dialog box.

The software can flag wells as the result of:

- Low allele frequency
- Small sample population
- Outliers mistakenly identified as a cluster (one or two samples in a cluster may in fact be an outlier)

## **Troubleshooting**

Review the calls and/or re-run the plate.

Notes		

6

## Thresholding Algorithm Failed (TAF)

## Description/ Function

This flag indicates that the software could not calculate a threshold because of a computational error.

**Note:** The Exponential Region, Baselining, Thesholding, and  $C_T$  Calculation Algorithms are performed in sequence. If one flag fails, then the SDS software does not compute the flags for the remaining algorithms in the sequence.

## **Troubleshooting**

- 1. Select the affected sample in the Results Table.
- **2.** Select the **Results** tab.
- **3.** In the Amplification Plot, review the amplification curves. If necessary, set the baseline and threshold manually for the affected run.

Notes			

# Adding Custom Dyes to the Pure Dye Set

The Applied Biosystems 7900HT Fast Real-Time PCR System can be used to run assays designed with custom dyes (dyes not manufactured by Applied Biosystems). However, before using custom dyes with the 7900HT Fast System, you must create and run a pure dye plate made with the custom dyes. The purpose of the custom pure dye plate is similar to that of a standard pure dye plate: the SDS software uses the custom pure dye plate to create a spectral standard to distinguish the custom dye.

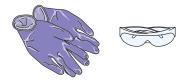
**IMPORTANT!** To use a custom dye on your 7900HT Fast System, it must fluoresce within the spectral range measured by the instrument. The working range of the CCD camera is 500 to 660 nm.

# **Before You Begin**

## Materials Required

You need the materials listed below to perform this procedure.

Powder-free gloves and safety goggles



Reaction plates appropriate for your sample block:

- Standard 384-well reaction plates
- Standard 96-well reaction plates
- Fast 96-well reaction plates

**Note:** You cannot create custom dyes for the TagMan<sup>®</sup> Low Density Array Block.

Pipettors and pipet tips (100- $\mu$ L)



# **Appendix A** *Before You Begin*

Water, deionized	
Custom dyes	
Tubes (2-mL and 10-mL)	
Optical adhesive covers or Optical Caps (PN 4323032)	
Centrifuge, with plate adapter	

## Workflow To add custom dyes to the pure dye set:

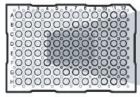
- **1.** Obtain the tools and materials required (see pages 141 and 142).
- **2.** Create a dilution series plate for each custom dye you want to add to the pure dye set (see page 144).
- **3.** Create a plate document for the dilution series plate (see page 144).
- **4.** Run and analyze the dilution series plate (see page 145).
- **5.** Create a pure dye plate with the custom dye(s) (see page 148).
- **6.** Add the custom dye(s) to the SDS software (see page 149).
- 7. Create a plate document template with the custom dye(s) (see page 150).

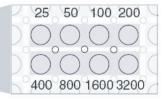
Notes			

# **Creating a Dilution Series Plate**

- 1. In the center wells of a reaction plate, prepare a dilution series of the custom dye (for example 25, 50, 100, 200, 400, 800, 1600, and 3200 nM concentration):
  - For a standard 384-well or Fast 96-well reaction plate, use 20 μL per well
  - For a standard 96-well reaction plate, use 50 μL per well

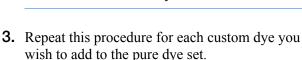






**2.** Seal the wells of the reaction plate using an optical adhesive cover or Optical Caps (PN 4323032).

**IMPORTANT!** Do not use MicroAmp® caps (domed) or Optical Tubes with the 7900HT Fast System. You can use Optical Caps (PN 4323032) (ONLY on the standard 96-well reaction plates) with the 7900HT Fast System.



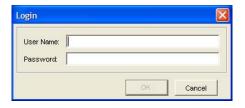


# Creating a Plate Document for the Dilution Series Plate

- Double-click on the desktop (or select Start > All Programs > Applied Biosystems > SDS 2.3 > SDS 2.3) to start the SDS software.
- 2. If the login option is enabled, the Login dialog box appears. Enter your User Name and Password, then click OK.

**Note:** If the login option is not enabled, no Login dialog box appears. Skip to step 3 below.

3. From the SDS software menu bar, click ☐ (or select File ➤ New).

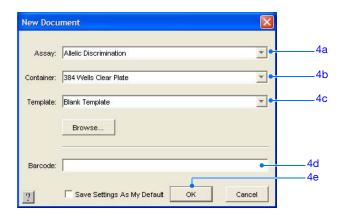


N.I	_	4	_	_
L/I	$\cap$	١T	$^{\Box}$	9

- **4.** Complete the New Document dialog box:
  - a. Assay Select Allelic Discrimination.
  - **b.** Container Select the appropriate format.
  - **c.** Template Select **Blank Template**.
  - d. Barcode Leave this field blank.
  - e. Click or The software creates and opens a new plate document.

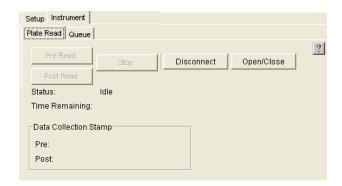
**Note:** It is not necessary to configure detector, sample, or method information for the dilution series plate document. The purpose of the run is to establish the correct working concentration for the dye by viewing the intensity of the raw spectra produced by the wells in the dilution series.

**5.** Repeat this procedure for each dilution series plate you created.



# Running and Analyzing the Dilution Series Plate

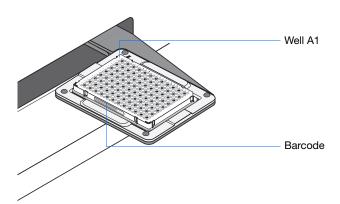
- 1. In the plate document, select the **Instrument** > **Plate Read** tabs.
- **2.** Click **Open/Close**. The instrument tray rotates to the OUT position.



**3.** Place the prepared series dilution plate into the instrument tray as shown.

**IMPORTANT!** The A1 position is located in the top-left side of the instrument.

#### Standard 384-well reaction plate



**4.** Click **Post Read.** The instrument tray rotates to the IN position and the instrument performs the run.

As the instrument performs the run, it displays status information in the Plate Read tab.

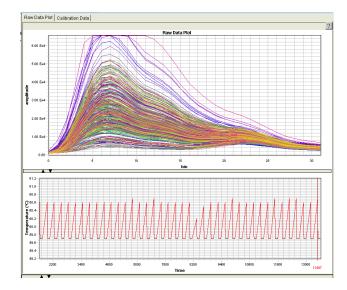
After the run, the status values and buttons are grayed-out, the Analysis button is enabled (), and a message indicates whether or not the run is successful.

- **5.** Click **Open/Close** to eject the reaction plate.
- **6.** Click (or select **Analysis** ▶ **Analyze**). The software analyzes the raw run data.
- 7. Click [3] (Show System Raw Data Pane).

**8.** In the Raw Data Plot, determine the highest concentration of dye that does not produce a saturated signal, then record it for future use.

Saturated signals are characterized by their high peaks that rise beyond detectable levels (> 65,000 fluorescent units) and appear as plateaus on the Raw Data plot.

The concentration of the custom dye that yields the highest possible signal but does not saturate is the maximum concentration for use with the 7900HT Fast System.



**9.** Repeat this procedure for each dilution series plate you created.

Notes			

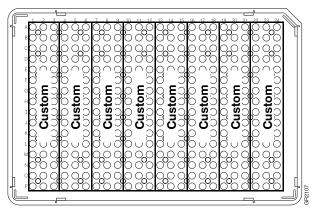
# Creating a Pure Dye Plate with the Custom Dye(s)

- **1.** Prepare 5 mL of a custom dye at the concentration determined in step 8 on page 147.
- **2.** Pipet the diluted custom dye to at least three columns of a reaction plate, as shown at right.

**IMPORTANT!** The optical configuration of the 7900HT instrument requires that each custom dye occupy at least three columns of the reaction plate to permit adequate data collection.

- For a standard 384-well or Fast 96-well reaction plate, pipet 20 µL per well
- For a standard 96-well reaction plate, pipet 50 µL per well





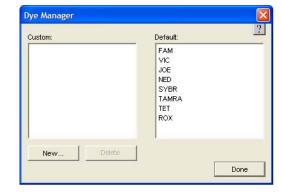
- **3.** Repeat steps 1 through 2 for each custom dye you wish to add to the reaction plate.
- **4.** Seal the wells of the reaction plate using an optical adhesive cover or Optical Caps (PN 4323032).

**IMPORTANT!** Do not use MicroAmp® caps (domed) or Optical Tubes with the 7900HT Fast System. You can use Optical Caps (PN 4323032) (ONLY on the standard 96-well reaction plates) with the 7900HT Fast System.



# Adding the New Custom Dye(s) to the SDS Software

- 1. From the SDS software menu bar, select **Tools** > **Dye Manager**.
- 2. In the Dye Manager dialog box, click New...



- **3.** In the Add Dye dialog box:
  - a. Enter a name for the custom dye.
  - b. Click ok The new dye appears in the Dye Manager's Custom dye list.

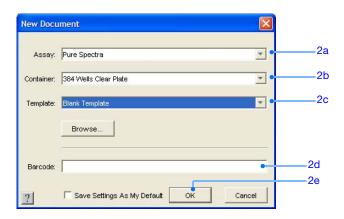


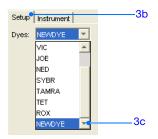


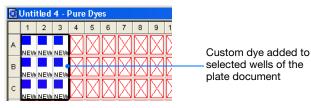
- **4.** Repeat steps 1 through 3 for each custom dye you added to the custom pure dye plate (page 148).
- **5.** Click Done The SDS software makes the new dye(s) available to pure dye plate documents.

# Creating a Plate Document Template with the Custom Dye(s)

- 1. Click □ (or select File > New).
- **2.** Complete the New Document dialog box:
  - a. Assay Select Pure Spectra.
  - **b.** Container Select the appropriate format.
  - c. Template Select Blank Template.
  - d. Barcode Leave this field blank.
  - e. Click \_\_\_ok\_\_\_. The software creates and opens a new plate document.
- **3.** Apply the new custom dye(s) to the plate document:
  - **a.** Select the wells containing the first custom dye.
  - **b.** Select the **Setup** tab.
  - **c.** In the Dyes drop-down list, select the appropriate dye. The software applies the dye to the selected wells.
  - **d.** Repeat steps a and b to configure the plate document with any additional custom dyes.



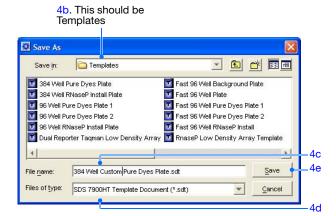




- **4.** Save the custom pure dye plate document as a plate document template:
  - a. Click (or select File > Save) to open the Save As dialog box.
  - b. In the Save in field, navigate to and open
     AppliedBiosystems ➤ SDS2.3 ➤
     Templates.

**Note:** By saving the plate document template to the Templates directory, it becomes available from the Template dropdown list in the New Document dialog box.

- **c.** In the File name field, enter a name for the plate document template.
- d. Select SDS 7900HT Template Document (\*.sdt) from the Files of type drop-down list.
- e. Click \_\_\_\_\_. The software saves the plate document as a plate document template.
- **5.** To run the custom dyes, see "Performing a Pure Dye Calibration" on page 17.



# Appendix A Creating a Plate Document Template with the Custom Dye(s) Notes

# **Instrument Connections**

**Electrical** Figure B-1 and Table B-1 illustrate the electrical connections of the Applied Biosystems 7900HT Fast Real-Time PCR System components.

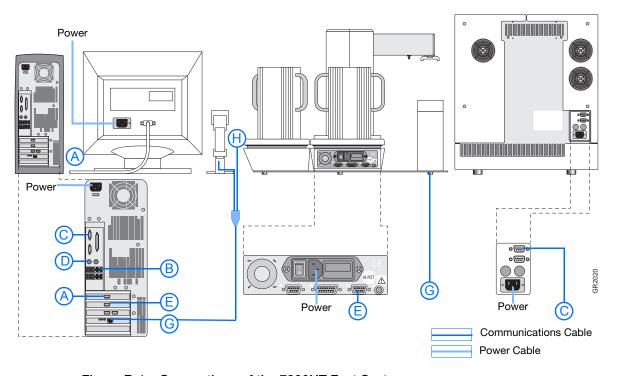


Figure B-1 Connections of the 7900HT Fast System

Table B-1 Connections of the 7900HT Fast System

Cable	Туре	Connects	То
Α	Communication	Computer (Monitor Port)	Monitor
В	Comm/Power	Computer (Mouse Port)	Mouse (not shown)
С	Serial	Computer (Serial Port 1)	7900HT Instrument
D	Comm/Power	Computer (Keyboard Port)	Hand-held Bar Code Reader
E	Communication	Computer (Serial Port 2)	Plate Handler (Port C)
G‡	Comm/Power	Computer (ISA Card 1)	Fixed-Position Bar Code Reader
Н	Comm/Power	Bar Code Reader Cable	Keyboard (not shown)

‡See Figure B-2 on page B-154.

## Fixed-Position Bar Code Reader

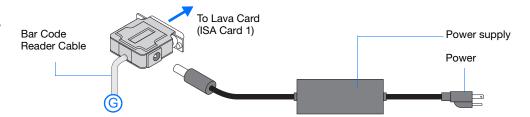


Figure B-2 Fixed-position bar code reader connection

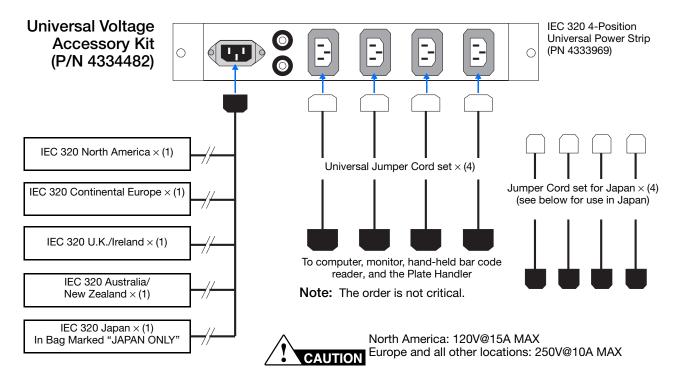


Figure B-3 Universal power strip

## To install the Universal Voltage Accessory Kit (All Countries Except Japan):

- **1.** Connect one universal jumper to each accessory (see below for use in Japan).
- **2.** Connect the other end of the power jumpers to the power strip outputs.
- **3.** Choose the correct country specific power input cord for the geographical region and connect it to the power strip input.
- **4.** Power off all instrument accessories (monitor, computer, and Plate Handler).
- **5.** Connect the power-input cord to the AC outlet.

- **6.** Power on accessories.
- **7.** Discard unused cord sets.

## Usage Guidelines for Japan

- **1.** Use only the cord sets supplied in the bag marked "JAPAN ONLY."
- **2.** Discard all other unused cord sets.

Notes			



Notes\_\_\_\_\_

# Parts and Consumables

**Note:** Part numbers listed in this appendix are for customers inside the United States. Contact your Regional Sales Office for local Part numbers and prices. See "How to Obtain Support" on page vii.

# Interchangeable Sample Block Modules and Accessories

The Applied Biosystems 7900HT Fast Real-Time PCR System features a Peltier-based, interchangeable sample block based on the technology established in the GeneAmp® PCR System 9700 thermal cycler.

The use of an interchangeable sample block:

- Reduces instrument downtime by allowing immediate replacement of the block.
- Permits easy access to the sample block for troubleshooting and maintenance.
- Supports multiple consumable formats.
- Provides several different modes of operation (including Max mode and programmable temperature ramps).

Table C-1 Sample block modules and accessories

Part No.	Description	Quantity
4331406	7900HT System Standard 384-Well Block Upgrade Kit	1 kit
	Includes a Standard 384-Well Block, a 384-Well Plate Adapter, and a Sequence Detection Systems 384-Well Spectral Calibration Kit (PN 4323977)	
4331405	7900HT System Standard 96-Well Block Upgrade Kit	1 kit
	Includes a Standard 96-Well Block, a 96-Well Plate Adapter, and an ABI PRISM® 7900HT Sequence Detection Systems 96-Well Spectral Calibration Kit (PN 4328639)	
4351402	7900HT System Fast 96-Well Block Upgrade Kit	1 kit
	Includes a Fast 96-Well Block, a Fast 96-Well Plate Adapter, a 7900HT System Fast 96-Well Spectral Calibration Kit (PN 4351653), a TaqMan® RNase P Fast 96-Well Instrument Verification Plate (PN 4351979), and a TaqMan® Fast Reagents Starter Kit (PN 4352407)	

Notes		

Table C-1 Sample block modules and accessories (continued)

Part No.	Description	Quantity
4329012	7900HT System TaqMan® Low Density Array Upgrade	1 kit
	Includes a TaqMan® Low Density Array Block, microfluidic card sealer, centrifuge buckets and adapters, and a chemistry installation kit.	

# **Consumables and Disposables**

The 7900HT Fast System can run:

- 384-Well Optical Reaction Plates (also referred to as *standard 384-well reaction plates*) sealed with optical adhesive covers
- MicroAmp® 96-Well Optical Reaction Plates (also referred to as standard 96-well reaction plates) sealed with optical adhesive covers or Optical Caps (PN 4323032, flat cap strips only)
- Optical 96-Well Fast Thermal Cycling Plates (also referred to as *Fast 96-well reaction plates*) sealed with optical adhesive covers
- TaqMan® Low Density Arrays

The optical plates recommended above are designed specifically for fluorescence-based PCR chemistries and are frosted to minimize external fluorescent contamination. Before running prepared optical plates on the 7900HT instrument, each plate must be sealed with the optical adhesive covers recommended above. Applied Biosystems optical adhesive covers are specifically designed to permit the transmission of light to and from the wells of the optical plate.

**IMPORTANT!** Do not use MicroAmp<sup>®</sup> caps (domed) or Optical Tubes with the 7900HT Fast System. You can use Optical Caps (PN 4323032) (ONLY on the standard 96-well reaction plates) with the 7900HT Fast System.

Table C-2 Consumables and disposables for the 7900HT instrument

Part No.	Description	Quantity
Optical Adh	nesive Covers	
4313663	ABI PRISM® Optical Adhesive Cover Starter Kit	20 Covers
	Includes 20 ABI PRISM® Optical Adhesive Covers, an Applicator, and an ABI PRISM® Optical Cover Compression Pad.	
4311971	ABI PRISM® Optical Adhesive Covers	100 Covers
4360954	Optical Adhesive Covers	25 Covers

11/1		

Table C-2 Consumables and disposables for the 7900HT instrument (continued)

Part No.	Description	Quantity
4323032	Optical Caps, 8 Caps/Strip	300 Strips/ Pkg
		2400 Caps/ Pkg
Standard 3	84-Well Reaction Plates	
4309849	384-Well Clear Optical Reaction Plate with Barcode (code 128)	50 Plates
4326270	384-Well Clear Optical Reaction Plate with Barcode (code 128), 10-Pack	500 Plates
	Includes 10 of PN 4309849, 384-Well Clear Optical Reaction Plates with Barcode	
Standard 9	6-Well Reaction Plates	
4306737	MicroAmp® 96-Well Optical Reaction Plate with Barcode (code 128)	20 Plates
4326659	MicroAmp® 96-Well Optical Reaction Plate with Barcode (code 128), 25-Pack	500 Plates
	Includes 25 of PN 4306737, MicroAmp® 96-Well Optical Reaction Plates with Barcode	
4314320	MicroAmp® 96-Well Optical Reaction Plate with Barcode (code 128) and ABI PRISM® Optical Adhesive Covers	100 Plates
	Includes 100 ABI PRISM® Optical Adhesive Covers (PN 4311971) and 5 of PN 4306737, MicroAmp® 96-Well Optical Reaction Plates with Barcode	
4312063	MicroAmp <sup>®</sup> Splash Free Support Base for 96-Well Reaction Plates	10 Bases
Fast 96-We	Il Reaction Plates	
4346906	Optical 96-Well Fast Thermal Cycling Plate with Barcode (code 128)	20 Plates
4312063	MicroAmp® Splash Free Support Base for 96-Well Reaction Plates	10 Bases
TaqMan® L	ow Density Arrays	
	TaqMan® Low Density Arrays	Variable*

<sup>\*</sup> See the Applied Biosystems Web site.

Notes\_\_\_\_

# **Instrument Maintenance and Verification**

The following sequence detection kits and reagents are used to perform routine maintenance on and verify the function of the 7900HT Fast System.

Table C-3 Consumables for calibration and verification runs

Part Number	Description	Quantity	
Sequence l	Detection Systems Calibration Kits		
4328639	ABI PRISM® 7900HT Sequence Detection Systems 96-Well Spectral Calibration Kit	3 x 96-Well Plates	
	Includes three MicroAmp <sup>®</sup> 96-Well Optical Reaction Plates: one preloaded and sealed background plate, and two preloaded and sealed pure dye plates containing eight separate dye standards (FAM <sup>™</sup> , JOE <sup>™</sup> , NED <sup>™</sup> , ROX <sup>™</sup> , SYBR <sup>®</sup> Green, TAMRA <sup>™</sup> , TET <sup>™</sup> , VIC <sup>®</sup> ).		
4323977	Sequence Detection Systems 384-Well Spectral Calibration Kit	2 x	
	Includes two 384-Well Optical Reaction Plates: one preloaded and sealed background plate and one preloaded and sealed pure dye plate containing eight separate dye standards (FAM, JOE, NED, ROX, SYBR Green, TAMRA, TET, VIC).	384-Well Plates	
4351653	7900HT System Fast 96-Well Spectral Calibration Kit	3 x 96-Well	
	Includes three Optical 96-Well Fast Thermal Cycling Plates: one preloaded and sealed background plate, and two preloaded and sealed pure dye plates containing eight separate dye standards (FAM™, JOE™, NED™, ROX™, SYBR® Green, TAMRA™, TET™, VIC®).	Plates	
4362745	TaqMan® Low Density Array Spectral Calibration Kit	4 x TLDAs	
	Includes four empty TaqMan <sup>®</sup> Low Density Arrays (TLDAs) and four tubes: one background solution, one FAM <sup>™</sup> dye, one ROX <sup>™</sup> dye, and one VIC <sup>®</sup> dye.		
Instrument	Verification Plates		
4310982	TaqMan® RNase P Instrument Verification Plate	1 x 96-Well	
	Includes one MicroAmp <sup>®</sup> 96-Well Optical Reaction Plate. Each well contains preloaded reaction mix (master mix, RNase P primers, and FAM <sup>™</sup> dye-labeled probe) and template to detect and quantitate genomic copies of the human RNase P gene.	Plate	
4323306	TaqMan® RNase P 384-Well Instrument Verification Plate	1 x 384-	
	Includes one 384-Well Optical Reaction Plate. Each well contains preloaded reaction mix (master mix, RNase P primers, and FAM <sup>™</sup> dye-labeled probe) and template to detect and quantitate genomic copies of the human RNase P gene.	Well Plate	

Notes			

Table C-3 Consumables for calibration and verification runs (continued)

Part Number	Description	Quantity
4351979	TaqMan <sup>®</sup> RNase P Fast 96-Well Instrument Verification Plate Includes one Optical 96-Well Fast Thermal Cycling Plate. Each well contains preloaded reaction mix (master mix, RNase P primers, and FAM <sup>™</sup> dye-labeled probe) and template to detect and quantitate genomic copies of the human RNase P gene.	1 x 96-Well Plate
4351468	TaqMan® Low Density Array RNase P Kit Includes one empty TaqMan® Low Density Array (TLDA) and eight tubes of solution. Each solution contains reaction mix (master mix, RNase P primers, and FAM™-MGB dye-labeled probe) and template to detect and quantitate genomic copies of the human RNase P gene.	1 TLDA

Appendix C Instrument Maintenance and Verification	
Notes	

# Index

A	bold text, when to use v
absolute quantification troubleshooting 124	С
adjusting plate-sensor switch 92 to 96 adjustment knob 89	calibrating the 7900HT instrument adjusting the plate-sensor switch 92 aligning the fixed-position bar code readers 110
air bubbles 121	aligning the Plate Handler 97
CLV fixed-position bar code reader 112 to 113 fixed-position bar code readers 110 LD fixed-position bar code reader 114 to 116 Plate Handler 97 to 108 amplification run starting 145 analyzing background data 14 pure dye data 29 Applied Biosystems contacting vii customer feedback on documentation vii Information Development department vii Technical Support vii assumptions for using this guide v Automation Accessory	calibration background 3, 17  CAUTION, description vi changing gripper finger pads 90 sample block module 53 chemistry non-optimized 121 troubleshooting 120 to 123 cleaning gripper finger pads 91 sample block wells 66, 67 computer troubleshooting 84 consumables Fast 96-well reaction plates 159 improper or damaged plastics 123 Optical Adhesive Covers 158
components 89 Plate Handler, See <i>Plate Handler</i> See <i>fixed-position bar code reader</i>	Optical Adhesive Covers 138 Optical Cap Strips 158 standard 384-well reaction plates 159 standard 96-well reaction plates 159 TaqMan Low Density Arrays 159 writing on reaction plates 122
background calibration about 3, 17 analyzing the data 14 creating the plate document 10 preparing the plate 6 running the plate 12 when to perform 3, 17	contamination decontaminating the sample block 65 fluorescent, common sources 122 isolating on the sample block module 15 conventions bold text v for describing menu commands v
background component, about 3 background plate creating 8 preparing 6	IMPORTANTS! vi in this guide v italic text v Notes v user attention words v
running 12 background run troubleshooting 15 backing up SDS files 73 backup storage devices 73	creating background plate 8 background plate document 10 custom pure dye plate documents 150 pure spectra plate document 24

verification run plate document 41	reader 112 to 113
custom pure dyes 141	launching 112, 114
customer feedback, on Applied Biosystems	LDHOST software
documents vii	aligning the LD fixed-position bar code reader 114 to 116
D	low copy templates 123
data management 73	M
decontaminating the sample block module 65 to 66	
documentation, related vi	maintenance schedule 52, 72, 88 calibration and verification runs 2
F	master mixes using 121
finger pads	menu commands, conventions for describing v
cleaning 91	modes of operation 74
replacing 90	automated 75
fixed-position bar code reader	stand-alone 74
aligning CLV 112 to 113 aligning LD 114 to 116	MSDSs, obtaining vi, vii
connections 153	P
fluorescent	performing
contamination 122 detection system 153	background calibration 6
detection system 133	pure dye run 21
G	pipetting errors 121
	pipettors, using 121
gripper 89	plate adapter, changing 62
	plate document
Н	background 10
hand-held bar code reader	pure spectra 24
connections 153	verification run 41
	Plate Handler 89
I control of the second of the	adjustment knob 89
imprecise pipetting 121	aligning 97 to 108 cleaning the finger pads 91
improper threshold setting 120	gripper 89
Information Development department, contacting vii	plate stack positions 89
installing	plate-sensor switch 89
plate adapter 62	replacing the finger pads 91
sample block module 53	plate stacks
SDS software 83	positions 89 plate-sensor switch 89
SDS Software Updates 83 Windows service pack updates 83	adjusting 92 to 96
instrument	precision
troubleshooting 84	causes of low precision 120 to 123
instrument tray	preparing
replacing the plate adapter 62	background plate 6
irreproducibility	pure dye plates 21
causes 120 to 123	RNase P plate 37
italic text, when to use v	RNase P TLDA 39
	process quality values (PQV)
L	rules for 129 pure dye calibration
LAVA software	analyzing the data 29
aligning the CLV fixed-position bar code	running the plate 27

pure dye plate	T
constructing for custom dyes 141	TaqMan Low Density Array RNase P Kit
running 27	about 34
pure dye run	TaqMan RNase P Instrument Verification Plates
troubleshooting 31	about 32
pure spectra calibration	analyzing 47
about 17	kits 160
creating the plate document 24	Technical Support, contacting vii
preparing pure dye plates 21	text conventions v
_	
R	threshold
reagents	improper setting 120
non-Applied Biosystems PCR reagents 123	TLDA RNase P card
TaqMan RNase P Instrument Verification	about 34
Plates 160	training, information on vii
reinstalling the SDS software 83	troubleshooting
relative quantification	7900HT instrument 84
troubleshooting 124	background runs 15
replacing	chemistry problems 120 to 123
gripper finger pads 90, 91	computer 84 end-point runs 126
sample block module 53	fixed-position bar code reader 117
RNase P plate	pure dye runs 31
running 43	real-time runs 124
RNase P TLDA	SDS software 84
running 43	Zymark Twister Microplate Handler 117
RNase P verification run	
preparing the plate 37	U
preparing the RNase P TLDA 39	undating
running	updating
background plate 12	SDS Software 83 Windows operating system 83
pure dye plate 27	user attention words, described v
RNase P plate 43	user attention words, described
RNase P TLDA 43	W
	V
S	verification run
comple blook leaking her 57	creating the plate document 41
sample block locking bar 57	running the RNase P plate or RNase P TLDA 43
sample block locking bolt 57	
sample block module	W
cleaning sample block module wells 66, 67	WARNING, description vi
contamination 122	-
replacing 53	Windows service pack updates, installing 83
saturation, signal 147	_
SDS software	Z
installing 83	Zymark Twister Software
reinstalling 83	aligning the Plate Handler 97 to 108
signal saturation 147	testing the plate sensor switch 95
software, starting 10, 24, 41, 56, 60, 63, 79, 81, 82, 99,	
111, 144	
spectral calibration, See pure dye runs	
storage devices, backup 73	

Index

#### Worldwide Sales and Support

Applied Biosystems vast distribution and service network, composed of highly trained support and applications personnel, reaches 150 countries on six continents. For sales office locations and technical support, please call our local office or refer to our Web site at www.appliedbiosystems.com.

Applied Biosystems is committed to providing the world's leading technology and information for life scientists.

#### Headquarters

850 Lincoln Centre Drive Foster City, CA 94404 USA Phone: +1 650.638.5800 Toll Free (In North America): +1 800.345.5224 Fax: +1 650.638.5884

06/2010

