

Applied Biosystems 7500/7500 Fast Real-Time PCR Systems System Maintenance



Applied Biosystems 7500/7500 Fast Real-Time PCR Systems System Maintenance

Overview

1

Perform the
Regions of Interest
(ROI) Calibration

2

Perform the
Background
Calibration and
Optical Calibration

3

Perform the Dye
Calibration

4

Verify the
Instrument
Performance

5

User-Performed
Maintenance

6

© Copyright 2008, 2010 Applied Biosystems. All rights reserved.

Information in this document is subject to change without notice. Applied Biosystems assumes no responsibility for any errors that may appear in this document.

APPLIED BIOSYSTEMS DISCLAIMS ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE. IN NO EVENT SHALL APPLIED BIOSYSTEMS BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

NOTICE TO PURCHASER: Label License

The Applied Biosystems 7500 and 7500 Fast Real-Time PCR Systems are 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:

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

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

SYBR is a registered trademark of Molecular Probes, Inc.

Microsoft and Windows are registered trademarks of Microsoft Corporation.

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

Part Number 4387777 Rev. D
06/2010

Contents

	Preface	vii
	How to Use This Guide	vii
	How to Obtain More Information	viii
	How to Obtain Support	x
	Safety and EMC Compliance Information	xi
	Safety Conventions Used in This Document	xii
	Symbols on Instruments	xiii
	Safety Labels on Instruments	xiv
	General Instrument Safety	xvi
	Chemical Safety	xvii
	Chemical Waste Safety	xix
	Electrical Safety	xx
	Physical Hazard Safety	xx
	Biological Hazard Safety	xxi
	Workstation Safety	xxi
	Safety and Electromagnetic Compatibility (EMC) Standards	xxii
Chapter 1	Overview	1
	About the 7500/7500 Fast System	2
	How to Use This Guide	4
	Recommended Maintenance Schedule	5
	Maintain the Computer Hard Drive(s)	6
	Archive and Back Up EDS Files	6
Chapter 2	Perform the Regions of Interest (ROI) Calibration	7
	Overview	8
	Prepare the ROI Calibration Plate	9
	Perform the Calibration	12
	Perform an Automated ROI Calibration	12
	Perform a Manual ROI Calibration	13
	Troubleshoot the ROI Calibration	17

Chapter 3	Perform the Background Calibration and Optical Calibration	19
	Perform the Background Calibration	20
	Prepare the Background Calibration Plate	21
	Perform the Background Calibration	23
	Perform the Optical Calibration	25
	Prepare the Calibration Plate	25
	Perform the Optical Calibration	28
	Troubleshoot the Background Calibration	30
Chapter 4	Perform the Dye Calibration	31
	Overview	32
	Prepare the Dye Calibration Plates	35
	Perform the Dye Calibration	36
	Troubleshoot the Dye Calibration	43
Chapter 5	Verify the Instrument Performance	45
	Overview	46
	Set Up the Experiment	48
	Run the Experiment	49
	Analyze the Experiment	51
	Troubleshoot the RNase P Experiment	54
Chapter 6	User-Performed Maintenance	55
	Monitor the 7500/7500 Fast System	56
	View the Instrument Log	56
	Monitor the Instrument Status	57
	Monitor the Lamp Status	58
	Decontaminate the Sample Block	59
	Replace the Halogen Lamp	63
	Replace the Instrument Fuses	66
	Update the Windows Operating System	67
	Update the 7500 Software	68
Appendix A	Store, Move, and Install the 7500/7500 Fast System . . .	69
	Store the 7500/7500 Fast System	70
	Move the 7500/7500 Fast System	71
	Set Up the 7500/7500 Fast System	73

Appendix B **Create a Custom Dye Plate** 75

Appendix C **Create a Background Plate** 81

Index 83

How to Use This Guide

Purpose of This Guide The *Applied Biosystems 7500/7500 Fast Real-Time PCR System Guide* provides the information you need to maintain your Applied Biosystems 7500/7500 Fast Real-Time PCR System. This manual is designed to supplement the:

- Applied Biosystems 7500/7500 Fast Real-Time PCR System Site Preparation Guide
- Applied Biosystems 7500/7500 Fast Real-Time PCR System Computer Setup Guide

Audience This guide is intended for novice and experienced 7500/7500 Fast system users who need to maintain their system.

Assumptions This guide assumes that your 7500/7500 Fast system has been installed by an Applied Biosystems technical representative and that you:

- Are familiar with the Microsoft® Windows® operating system.
- Understand general techniques for preparing and handling DNA samples.
- Have a general understanding of hard drives and data storage, file transfers, and copying and pasting.

Text Conventions This guide uses the following conventions:

- **Bold** text 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 symbol (▶) separates successive commands you select from a drop-down or shortcut menu. For example:
Select **File ▶ Open ▶ Spot Set**.
Right-click the sample row, then select **View Filter ▶ View All Runs**.

User Attention Words Two 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.

Examples of the user attention words appear below:

Note: The Calibrate function is also available in the Control Console.

IMPORTANT! To verify your client connection to the database, you need a valid user ID and password.

Safety Alert Words Safety alert words also appear in user documentation. For more information, see “[Safety Alert Words](#)” on page xii.

How to Obtain More Information

Related Documentation The following related documents are shipped with the 7500/7500 Fast system:

Guide	Purpose and Audience	PN
<i>Applied Biosystems 7500/7500 Fast Real-Time PCR System Getting Started Guide for Genotyping Experiments</i>	Explains how to perform experiments on the 7500/7500 Fast system. Each Getting Started Guide functions as both: <ul style="list-style-type: none"> • Tutorial, using example experiment data provided with the Applied Biosystems 7500/7500 Fast Real-Time PCR Software (7500 Software). • Guide for your own experiments. Intended for laboratory staff and principal investigators who perform experiments using the 7500/7500 Fast system.	4387784
<i>Applied Biosystems 7500/7500 Fast Real-Time PCR System Getting Started Guide for Presence/Absence Experiments</i>		4387785
<i>Applied Biosystems 7500/7500 Fast Real-Time PCR System Getting Started Guide for Relative Standard Curve and Comparative C_T Experiments</i>		4387783
<i>Applied Biosystems 7500/7500 Fast Real-Time PCR System Getting Started Guide for Standard Curve Experiments</i>		4387779
<i>Applied Biosystems 7500/7500 Fast Real-Time PCR System Maintenance Guide</i>	Explains how to install and maintain the 7500/7500 Fast system. Intended for laboratory staff responsible for the installation and maintenance of the 7500/7500 Fast system.	4387777
<i>Applied Biosystems 7500/7500 Fast Real-Time PCR System Computer Setup Guide</i>		4387778
<i>Applied Biosystems 7500/7500 Fast Real-Time PCR System Reagent Guide</i>	Provides information about the reagents you can use on the 7500/7500 Fast system, including: <ul style="list-style-type: none"> • An introduction to TaqMan® and SYBR® Green reagents • Descriptions and design guidelines for the following experiment types: <ul style="list-style-type: none"> – Quantitation experiments – Genotyping experiments – Presence/absence experiments Intended for laboratory staff and principal investigators who perform experiments using the 7500/7500 Fast system.	4387787

Guide	Purpose and Audience	PN
<i>Applied Biosystems 7500/7500 Fast Real-Time PCR System Site Preparation Guide</i>	Explains how to prepare your site to receive and install the 7500/7500 Fast system. Intended for personnel who schedule, manage, and perform the tasks required to prepare your site for installation of the 7500/7500 Fast system.	4387776
<i>Applied Biosystems 7500/7500 Fast Real-Time PCR Software v2.0 Help</i>	Explains how to use the 7500 Software to: <ul style="list-style-type: none"> • Set up, run, and analyze experiments using the 7500/7500 Fast system. • Monitor a networked 7500/7500 Fast instrument. • Calibrate a 7500/7500 Fast instrument. • Verify the performance of a 7500/7500 Fast instrument with an RNase P run. Intended for: <ul style="list-style-type: none"> • Laboratory staff and principal investigators who perform experiments using the 7500/7500 Fast system. • Laboratory staff responsible for the installation and maintenance of the 7500/7500 Fast system. 	NA

Note: To open the user documentation included on the Documentation CD, use the Adobe® Acrobat® Reader® software available from www.adobe.com.

Note: For additional documentation, see “How to Obtain Support” on page x.

Obtaining Information from the Help System

The 7500 software has a Help system that describes how to use each feature of the user interface. Access the Help system by doing one of the following:

- Click  in the toolbar of the 7500 software window
- Select **Help ▶ 7500 Software Help**
- Press **F1**

You can use the Help system to find topics of interest by:

- Reviewing the table of contents
- Searching for a specific topic
- Searching an alphabetized index

You can also access PDF versions of all documents in the Applied Biosystems 7500/7500 Fast Real-Time PCR System document set from the Help system.

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

IMPORTANT! The e-mail address above is only for submitting comments and suggestions relating to documentation. To order documents, download PDF files, or for help with a technical question, go to www.appliedbiosystems.com, then click the link for **Support**. (See “How to Obtain Support” below).

How to Obtain Support

For the latest services and support information for all locations, go to www.appliedbiosystems.com, then click the link for **Support**.

At the Support page, you can:

- 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

In addition, the Support page provides access to worldwide telephone and fax numbers to contact Applied Biosystems Technical Support and Sales facilities.

Safety and EMC Compliance Information

This section includes the following topics:

■ Safety Conventions Used in This Document	xii
■ Symbols on Instruments	xiii
■ Safety Labels on Instruments	xiv
■ General Instrument Safety	xvi
■ Chemical Safety	xvii
■ Chemical Waste Safety	xix
■ Electrical Safety	xx
■ Physical Hazard Safety	xx
■ Biological Hazard Safety	xxi
■ Workstation Safety	xxi
■ Safety and Electromagnetic Compatibility (EMC) Standards	xxii

Safety Conventions Used in This Document

Safety Alert Words

Four safety alert words appear in Applied Biosystems user documentation at points in the document where you need to be aware of relevant hazards. Each alert word—**IMPORTANT**, **CAUTION**, **WARNING**, **DANGER**—implies a particular level of observation or action, as defined below:

Definitions

IMPORTANT! – Indicates 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.

 **DANGER** – Indicates an imminently hazardous situation that, if not avoided, will result in death or serious injury. This signal word is to be limited to the most extreme situations.

Except for **IMPORTANT**s, each safety alert word in an Applied Biosystems document appears with an open triangle figure that contains a hazard symbol. *These hazard symbols are identical to the hazard icons that are affixed to Applied Biosystems instruments (see “Safety Symbols” on page xiii).*

Examples

The following examples show the use of safety alert words:

IMPORTANT! Wear powder-free gloves when you handle the halogen lamp.

 **CAUTION** The lamp is extremely hot. Do not touch the lamp until it has cooled to room temperature.

 **WARNING** **CHEMICAL HAZARD.** Ethanol is a flammable liquid and vapor. Exposure causes eye, skin, and respiratory tract irritation and may cause central nervous system depression and liver damage. Read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

 **DANGER** **ELECTRICAL HAZARD.** Failure to ground the instrument properly can lead to an electrical shock. Ground the instrument according to the provided instructions.

Symbols on Instruments

Electrical Symbols on Instruments

The following table describes the electrical symbols that may be displayed on Applied Biosystems instruments.

Symbol	Description	Symbol	Description
	Indicates the On position of the main power switch.		Indicates a terminal that may be connected to the signal ground reference of another instrument. This is not a protected ground terminal.
	Indicates the Off position of the main power switch.		Indicates a protective grounding terminal that must be connected to earth ground before any other electrical connections are made to the instrument.
	Indicates a standby switch by which the instrument is switched on to the Standby condition. Hazardous voltage may be present if this switch is on standby.		Indicates a terminal that can receive or supply alternating current or voltage.
	Indicates the On/Off position of a push-push main power switch.		Indicates a terminal that can receive or supply alternating or direct current or voltage.

Safety Symbols

The following table describes the safety symbols that may be displayed on Applied Biosystems instruments. Each symbol may appear by itself or in combination with text that explains the relevant hazard (see [“Safety Labels on Instruments”](#) on [page xiv](#)). These safety symbols may also appear next to DANGERS, WARNINGS, and CAUTIONS that occur in the text of this and other product-support documents.

Symbol	Description
	Indicates that you should consult the manual for further information and to proceed with appropriate caution.
	Indicates the presence of an electrical shock hazard and to proceed with appropriate caution.
	Indicates the presence of a hot surface or other high-temperature hazard and to proceed with appropriate caution.
	Indicates the presence of a laser inside the instrument and to proceed with appropriate caution.
	Indicates the presence of moving parts and to proceed with appropriate caution.

Environmental Symbols on Instruments

The following symbol applies to all Applied Biosystems electrical and electronic products placed on the European market after August 13, 2005.

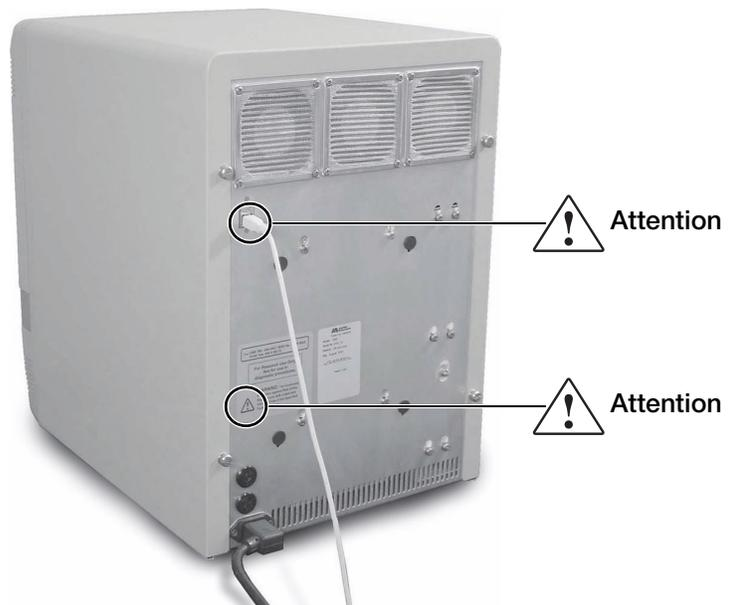
Symbol	Description
	<p>Do not dispose of this product as unsorted municipal waste. Follow local municipal waste ordinances for proper disposal provisions to reduce the environmental impact of waste electrical and electronic equipment (WEEE).</p> <p>European Union customers: Call your local Applied Biosystems Customer Service office for equipment pick-up and recycling. See http://www.appliedbiosystems.com for a list of customer service offices in the European Union.</p>

Safety Labels on Instruments

The following CAUTION, WARNING, and DANGER statements may be displayed on Applied Biosystems instruments in combination with the safety symbols described in the preceding section.

English	Français
CAUTION Hazardous chemicals. Read the Material Safety Data Sheets (MSDSs) before handling.	ATTENTION Produits chimiques dangereux. Lire les fiches techniques de sûreté de matériels avant la manipulation des produits.
CAUTION Hazardous waste. Refer to MSDS(s) and local regulations for handling and disposal.	ATTENTION Déchets dangereux. Lire les fiches techniques de sûreté de matériels et la réglementation locale associées à la manipulation et l'élimination des déchets.
WARNING Hot lamp.	AVERTISSEMENT Lampe brûlante.
WARNING Hot. Replace lamp with an Applied Biosystems lamp.	AVERTISSEMENT Composants brûlants. Remplacer la lampe par une lampe Applied Biosystems.
CAUTION Hot surface.	ATTENTION Surface brûlante.
DANGER High voltage.	DANGER Haute tension.
WARNING To reduce the chance of electrical shock, do not remove covers that require tool access. No user-serviceable parts are inside. Refer servicing to Applied Biosystems qualified service personnel.	AVERTISSEMENT Pour éviter les risques d'électrocution, ne pas retirer les capots dont l'ouverture nécessite l'utilisation d'outils. L'instrument ne contient aucune pièce réparable par l'utilisateur. Toute intervention doit être effectuée par le personnel de service qualifié de Applied Biosystems.
CAUTION Moving parts.	ATTENTION Parties mobiles.
WARNING This instrument is designed for 12V, 75W Halogen lamps only.	AVERTISSEMENT Cet instrument est conçu pour des lampes d'halogène de 12V et 75W seulement.

Locations of Warnings The Applied Biosystems 7500/7500 Fast Real-Time PCR System contains warnings at the locations shown below.



General Instrument Safety

 **WARNING** **PHYSICAL INJURY HAZARD.** Use this product only as specified in this document. Using this instrument in a manner not specified by Applied Biosystems may result in personal injury or damage to the instrument.

Moving and Lifting the Instrument

 **CAUTION** **PHYSICAL INJURY HAZARD.** The instrument is to be moved and positioned only by the personnel or vendor specified in the applicable site preparation guide. If you decide to lift or move the instrument after it has been installed, do not attempt to lift or move the instrument without the assistance of others, the use of appropriate moving equipment, and proper lifting techniques. Improper lifting can cause painful and permanent back injury. Depending on the weight, moving or lifting an instrument may require two or more persons.

Moving and Lifting Stand-Alone Computers and Monitors

 **WARNING** Do not attempt to lift or move the computer or the monitor without the assistance of others. Depending on the weight of the computer and/or the monitor, moving them may require two or more people.

Things to consider before lifting the computer and/or the monitor:

- Make sure that you have a secure, comfortable grip on the computer or the monitor when lifting.
- Make sure that the path from where the object is to where it is being moved is clear of obstructions.
- Do not lift an object and twist your torso at the same time.
- Keep your spine in a good neutral position while lifting with your legs.
- Participants should coordinate lift and move intentions with each other before lifting and carrying.
- Instead of lifting the object from the packing box, carefully tilt the box on its side and hold it stationary while someone slides the contents out of the box.

Operating the Instrument

Ensure that anyone who operates the instrument has:

- Received instructions in both general safety practices for laboratories and specific safety practices for the instrument.
- Read and understood all applicable Material Safety Data Sheets (MSDSs). See [“About MSDSs” on page xvii](#).

 **WARNING** **PHYSICAL INJURY HAZARD.** Use this instrument as specified by Applied Biosystems. Using this instrument in a manner not specified by Applied Biosystems may result in personal injury or damage to the instrument.

Cleaning or Decontaminating the Instrument

 **CAUTION** Before using a cleaning or decontamination method other than those recommended by the manufacturer, verify with the manufacturer that the proposed method will not damage the equipment.

Chemical Safety

Chemical Hazard Warning



WARNING CHEMICAL HAZARD. Before handling any chemicals, refer to the Material Safety Data Sheet (MSDS) provided by the manufacturer, and observe all relevant precautions.



WARNING CHEMICAL HAZARD. All chemicals in the instrument, including liquid in the lines, are potentially hazardous. Always determine what chemicals have been used in the instrument before changing reagents or instrument components. Wear appropriate eyewear, protective clothing, and gloves when working on the instrument.



WARNING CHEMICAL STORAGE HAZARD. Never collect or store waste in a glass container because of the risk of breaking or shattering. Reagent and waste bottles can crack and leak. Each waste bottle should be secured in a low-density polyethylene safety container with the cover fastened and the handles locked in the upright position. Wear appropriate eyewear, clothing, and gloves when handling reagent and waste bottles.

About MSDSs

Chemical manufacturers supply current Material Safety Data Sheets (MSDSs) with shipments of hazardous chemicals to *new* customers. They also provide MSDSs with the first shipment of a hazardous chemical to a customer after an MSDS has been updated. MSDSs provide the safety information you need to store, handle, transport, and dispose of the chemicals safely.

Each time you receive a new MSDS packaged with a hazardous chemical, be sure to replace the appropriate MSDS in your files.

Obtaining MSDSs

The MSDS for any chemical supplied by Applied Biosystems is available to you 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

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

Chemical Safety Guidelines

To minimize the hazards of chemicals:

- Read and understand the Material Safety Data Sheets (MSDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. (See “About MSDSs” on page xvii.)
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing). For additional safety guidelines, consult the MSDS.
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood). For additional safety guidelines, consult the MSDS.
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer’s cleanup procedures as recommended on the MSDS.
- Comply with all local, state/provincial, or national laws and regulations related to chemical storage, handling, and disposal.

Chemical Waste Safety

Chemical Waste Hazard



CAUTION HAZARDOUS WASTE. Refer to Material Safety Data Sheets and local regulations for handling and disposal.



WARNING CHEMICAL WASTE HAZARD. Wastes produced by Applied Biosystems instruments are potentially hazardous and can cause injury, illness, or death.



WARNING CHEMICAL STORAGE HAZARD. Never collect or store waste in a glass container because of the risk of breaking or shattering. Reagent and waste bottles can crack and leak. Each waste bottle should be secured in a low-density polyethylene safety container with the cover fastened and the handles locked in the upright position. Wear appropriate eyewear, clothing, and gloves when handling reagent and waste bottles.

Chemical Waste Safety Guidelines

To minimize the hazards of chemical waste:

- Read and understand the Material Safety Data Sheets (MSDSs) provided by the manufacturers of the chemicals in the waste container before you store, handle, or dispose of chemical waste.
- Provide primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary container. Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing). For additional safety guidelines, consult the MSDS.
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood). For additional safety guidelines, consult the MSDS.
- Handle chemical wastes in a fume hood.
- After emptying the waste container, seal it with the cap provided.
- Dispose of the contents of the waste tray and waste bottle in accordance with good laboratory practices and local, state/provincial, or national environmental and health regulations.

Waste Disposal

If potentially hazardous waste is generated when you operate the instrument, you must:

- Characterize (by analysis if necessary) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure the health and safety of all personnel in your laboratory.
- Ensure that the instrument waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.

IMPORTANT! Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.

Electrical Safety

 **DANGER ELECTRICAL SHOCK HAZARD.** Severe electrical shock can result from operating the Applied Biosystems 7500/7500 Fast Real-Time PCR System without its instrument panels in place. Do not remove instrument panels. High-voltage contacts are exposed when instrument panels are removed from the instrument.

Fuses  **WARNING FIRE HAZARD.** Improper fuses or high-voltage supply can damage the instrument wiring system and cause a fire. Before turning on the instrument, verify that the fuses are properly installed and that the instrument voltage matches the power supply in your laboratory.

 **WARNING FIRE HAZARD.** For continued protection against the risk of fire, replace fuses only with fuses of the type and rating specified for the instrument.

Power  **DANGER ELECTRICAL HAZARD.** Grounding circuit continuity is vital for the safe operation of equipment. Never operate equipment with the grounding conductor disconnected.

 **DANGER ELECTRICAL HAZARD.** Use properly configured and approved line cords for the voltage supply in your facility.

 **DANGER ELECTRICAL HAZARD.** Plug the system into a properly grounded receptacle with adequate current capacity.

Overvoltage Rating The Applied Biosystems 7500/7500 Fast Real-Time PCR System has an installation (overvoltage) category of II, and is classified as portable equipment.

Physical Hazard Safety

Moving Parts  **WARNING PHYSICAL INJURY HAZARD.** Moving parts can crush and cut. Keep hands clear of moving parts while operating the instrument. Disconnect power before servicing the instrument.

Biological Hazard Safety

General Biohazard



WARNING BIOHAZARD. Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. Follow all applicable local, state/provincial, and/or national regulations. Wear appropriate protective equipment, which includes but is not limited to: protective eyewear, face shield, clothing/lab coat, and gloves. All work should be conducted in properly equipped facilities using the appropriate safety equipment (for example, physical containment devices). Individuals should be trained according to applicable regulatory and company/institution requirements before working with potentially infectious materials. Read and follow the applicable guidelines and/or regulatory requirements in the following:

- U.S. Department of Health and Human Services guidelines published in *Biosafety in Microbiological and Biomedical Laboratories* (stock no. 017-040-00547-4; bml.od.nih.gov)
- Occupational Safety and Health Standards, Bloodborne Pathogens (29 CFR §1910.1030; www.access.gpo.gov/nara/cfr/waisidx_01/29cfr1910a_01.html).
- Your company's/institution's Biosafety Program protocols for working with/handling potentially infectious materials.

Additional information about biohazard guidelines is available at:

www.cdc.gov

Workstation Safety

Correct ergonomic configuration of your workstation can reduce or prevent effects such as fatigue, pain, and strain. Minimize or eliminate these effects by configuring your workstation to promote neutral or relaxed working positions.



CAUTION MUSCULOSKELETAL AND REPETITIVE MOTION HAZARD. These hazards are caused by potential risk factors that include but are not limited to repetitive motion, awkward posture, forceful exertion, holding static unhealthy positions, contact pressure, and other workstation environmental factors.

To minimize musculoskeletal and repetitive motion risks:

- Use equipment that comfortably supports you in neutral working positions and allows adequate accessibility to the keyboard, monitor, and mouse.
- Position the keyboard, mouse, and monitor to promote relaxed body and head postures.

Safety and Electromagnetic Compatibility (EMC) Standards

This section provides information on:

- U.S. and Canadian Safety Standards
- Canadian EMC Standard
- European Safety and EMC Standards
- Australian EMC Standards

U.S. and Canadian Safety Standards



This instrument has been tested to and complies with standard UL 61010A-1, “Safety Requirements for Electrical Equipment for Laboratory Use, Part 1: General Requirements” and with standard UL 61010-2-010, “Particular Requirements for Laboratory Equipment for the Heating of Materials.”

This instrument has been tested to and complies with standard CSA 1010.1, “Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use, Part 1: General Requirements.”

Canadian EMC Standard

This instrument has been tested to and complies with ICES-001, Issue 3: Industrial, Scientific, and Medical Radio Frequency Generators.

European Safety and EMC Standards



Safety

This instrument meets European requirements for safety (Low Voltage Directive 2006/95/EC). This instrument has been tested to and complies with standards EN 61010-1:2001, “Safety Requirements for Electrical Equipment for Measurement, Control and Laboratory Use, Part 1: General Requirements” and EN 61010-2-010, “Particular Requirements for Laboratory Equipment for the Heating of Materials,” and with standard EN 61010-2-081:2002+A1:2003 “Particular Requirements for Automatic and Semi-Automatic Laboratory Equipment for Analysis and Other Purposes.”

EMC

This instrument meets European requirements for emission and immunity (EMC Directive 2004/108/EC). This instrument has been tested to and complies with standard EN 61326 (Group 1, Class B), “Electrical Equipment for Measurement, Control and Laboratory Use – EMC Requirements.”

Australian EMC Standards



This instrument has been tested to and complies with standard AS/NZS 2064, “Limits and Methods Measurement of Electromagnetic Disturbance Characteristics of Industrial, Scientific, and Medical (ISM) Radio-frequency Equipment.”



Overview

1

This chapter covers:

- About the 7500/7500 Fast System 2
- How to Use This Guide 4
- Recommended Maintenance Schedule 5
- Maintain the Computer Hard Drive(s) 6

Note: For more information about any of the topics discussed in this guide, access the Help from within Applied Biosystems 7500/7500 Fast Real-Time PCR System Software v2.0 by pressing **F1**, clicking  in the toolbar, or selecting **Help ▶ 7500 Software Help**.

Notes

About the 7500/7500 Fast System

The Applied Biosystems 7500/7500 Fast Real-Time PCR System uses fluorescence-based polymerase chain reaction (PCR) reagents to provide:

- Quantitative detection of target nucleic acid sequences (targets) using real-time analysis.
- Qualitative detection of targets using post-PCR (endpoint) analysis.
- Qualitative analysis of the PCR product (achieved by melt-curve analysis that occurs post-PCR).

About Data Collection

The 7500/7500 Fast system collects raw fluorescence data at different points during a PCR, depending on the type of run that the instrument performs:

Run Type		Data Collection Point
Real-time runs	Standard curve	The instrument collects data following each extension step of the PCR.
	Relative standard curve	
	Comparative C _T ($\Delta\Delta C_T$)	
Post-PCR (endpoint) runs	Genotyping	The instrument collects data: <ul style="list-style-type: none"> • Before the PCR (For presence/absence experiments, data collection before the PCR is optional but recommended.) • (Optional) During the PCR. The instrument can collect data during the run (real-time); collecting data during the run can be helpful for troubleshooting. • After the PCR.
	Presence/Absence	

Regardless of the run type, a data collection point or *read* on the 7500/7500 Fast instrument consists of three phases:

- 1. Excitation** – The instrument illuminates all wells of the reaction plate, exciting the fluorophores in each reaction.
- 2. Emission** – The instrument optics collect the residual fluorescence emitted from the wells of the reaction plate. The resulting image consists only of light that corresponds to the range of emission wavelengths.
- 3. Collection** – The instrument assembles a digital representation of the residual fluorescence collected over a fixed time interval, then stores the raw fluorescence image for analysis.

After a run, the 7500 software uses regions of interest (ROI), optical, dye, and background calibration data to determine the location and intensity of the fluorescence signals in each read, the dye associated with each fluorescence signal, and the significance of the signal.

Notes _____

About the Filters The 7500/7500 Fast system uses five filters to support the following system dyes:

Filter 1	Filter 2	Filter 3	Filter 4	Filter 5
<ul style="list-style-type: none"> FAM™ dye SYBR® Green dye 	<ul style="list-style-type: none"> JOE™ dye VIC® dye 	<ul style="list-style-type: none"> TAMRA™ dye NED™ dye CY3® dye 	<ul style="list-style-type: none"> ROX™ dye Texas Red® dye 	Cy5® dye

For More Information

For information on

- The 7500/7500 Fast system – Refer to the *Applied Biosystems 7500/7500 Fast Real-Time PCR System Software Help*.

Note: To access the Help, select **Help ▶ 7500 Software Help** from within the 7500 software.

- Genotyping experiments – Refer to *Applied Biosystems 7500/7500 Fast Real-Time PCR System Getting Started Guide for Genotyping Experiments*.
- Presence/absence experiments – Refer to *Applied Biosystems 7500/7500 Fast Real-Time PCR System Getting Started Guide for Presence/Absence Experiments*.
- Relative standard curve and/or comparative C_T (C_T) experiments – Refer to *Applied Biosystems 7500/7500 Fast Real-Time PCR System Getting Started Guide for Relative Standard Curve and Comparative C_T Experiments*.
- Standard curve experiments – Refer to *Applied Biosystems 7500/7500 Fast Real-Time PCR System Getting Started Guide for Standard Curve Experiments*.

Notes _____

How to Use This Guide

This guide describes how to maintain the Applied Biosystems 7500/7500 Fast Real-Time PCR System. Chapters 2 through 5 of this manual describe calibrations that you must perform as regular maintenance of the 7500/7500 Fast system. Chapter 6 and the appendices contain maintenance procedures that you may need to resolve infrequent problems.

Chapter/ Appendix	Title	Description
2	Perform the Regions of Interest (ROI) Calibration	Describes how to perform an ROI calibration, which allows the 7500 software to map the positions of the wells on the sample block so that it can associate the fluorescence collected during a run with specific wells of the plate.
3	Perform the Background Calibration and Optical Calibration	Describes how to perform background and optical calibrations where the: <ul style="list-style-type: none"> Background calibration allows the 7500 software to remove the background fluorescence from experiment data. Optical calibration compensates for the physical effects of the fifth filter present in 7500/7500 Fast systems.
4	Perform the Dye Calibration	Describes how to perform dye calibrations, which allow the software to distinguish the individual contribution of each dye in the total fluorescence collected by the instrument.
5	Verify the Instrument Performance	Describes how to perform a TaqMan® RNase P Instrument Verification Plate run that can be used to verify the performance of a 7500/7500 Fast system.
6	User-Performed Maintenance	Describes how to: <ul style="list-style-type: none"> Replace the user-serviceable parts of the 7500/7500 Fast system. Resolve infrequent problems that can occur during instrument use.
A	Store, Move, and Install the 7500/7500 Fast System	Describes how to store, move, and reinstall the components of the 7500/7500 Fast system.
B	Create a Custom Dye Plate	Describes how to create a dye plate that can be used to calibrate the 7500/7500 Fast system for a dye not manufactured by Applied Biosystems.
C	Create a Background Plate	Describes how to create a background plate in the event that one is unavailable.

Notes _____

Recommended Maintenance Schedule

The following table displays the recommended maintenance schedule for the 7500/7500 Fast instrument and computer. The procedures listed in the table are intended for the user(s) of the 7500/7500 Fast system. To ensure proper operation of your instrument, perform the regular weekly, monthly, and semiannual maintenance indicated below.

IMPORTANT! The numbered lists in the table below indicate that the tasks must be performed in sequence.

Perform...	User-Performed Maintenance Task	See Page
Weekly	<ul style="list-style-type: none"> Check the computer disk space. If necessary, archive or back up your experiment files. Power off the computer controlling the 7500/7500 Fast instrument, then after 30 sec, power on the computer. Clean the surface of the 7500/7500 Fast instrument with a lint-free cloth. <p>IMPORTANT! Do not use organic solvents to clean the 7500/7500 Fast system.</p>	<p>6</p> <p>—</p> <p>—</p>
Monthly	<ol style="list-style-type: none"> Check the lamp status. If necessary, replace the halogen lamp. Perform a background calibration.† Run disk cleanup and disk defragmentation. 	<p>58</p> <p>20</p> <p>6</p>
Semiannually (6 Months)	<ol style="list-style-type: none"> Check the lamp status. If necessary, replace the halogen lamp. Perform a regions of interest (ROI) calibration. Perform a background calibration. Perform an optical calibration. Perform a dye calibration. Perform an RNase P instrument verification run. 	<p>58</p> <p>7</p> <p>20</p> <p>25</p> <p>31</p> <p>45</p>
As needed	<ul style="list-style-type: none"> Decontaminate the 7500/7500 Fast instrument. Replace the Halogen Lamp. Replace the 7500/7500 Fast instrument fuses. Update the Windows operating system. Update the 7500 software. 	<p>59</p> <p>63</p> <p>66</p> <p>67</p> <p>68</p>

† You can perform a background calibration to check for contamination. If any parts of the optics are replaced or moved, you must run an ROI calibration, a background calibration, an optical calibration, a dye calibration, and an RNase P instrument verification run.

Notes

Maintain the Computer Hard Drive(s)

When to Defragment and Clean Up the Hard Drive

- At least once every month
- When a message is displayed by the Windows operating system instructing you to defragment

For More Information

In the desktop, select **Start ▶ Help and Support** to access the Help for the Windows operating system. Use the search function of the Help to find information on the “Disk Cleanup” and “Disk Defragment” utilities.

IMPORTANT! Do not run the disk management utilities and 7500 software at the same time.

Archive and Back Up EDS Files

Archive EDS Files Regularly

To conserve space on the computer hard drive, older EDS files can be archived using a data compression utility. Several commercially available compression utilities are available. PKZIP and *.arc are archive formats common to the Microsoft® Windows® operating system.

Back Up EDS Files

Applied Biosystems strongly recommends that you back up your experiments.

Backing up data:

- Protects against potential loss of data caused by an unforeseen failure of the computer or its hard drive(s).
- Conserves space on the hard drive and optimizes performance, if you remove old data after backing up.

Develop a Data Management Strategy

Applied Biosystems recommends developing a strategy for managing the files produced by the 7500 software.

Note: Real-time runs generate significantly more data than genotyping or presence/absence experiments. During one day of real-time operation, the 7500/7500 Fast system can generate more than 10 MB of data.

Check Disk Space

If you perform real-time experiments on your 7500/7500 Fast system, check the amount of available space on your hard drive weekly. When the hard drive is within 20% of maximum capacity, transfer the older data to a backup storage device.

Notes _____

2

Perform the Regions of Interest (ROI) Calibration

This chapter covers:

■ Overview.....	8
■ Prepare the ROI Calibration Plate.....	9
■ Perform the Calibration	12
■ Perform an Automated ROI Calibration.....	12
■ Perform a Manual ROI Calibration	13
■ Troubleshoot the ROI Calibration.....	17

Note: For more information about any of the topics discussed in this guide, access the Help from within Applied Biosystems 7500/7500 Fast Real-Time PCR System Software v2.0 by pressing **F1**, clicking  in the toolbar, or selecting **Help ▶ 7500 Software Help**.

Notes

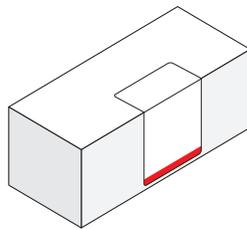
Overview

A regions of interest (ROI) calibration maps the positions of the wells on the sample block of the Applied Biosystems 7500/7500 Fast Real-Time PCR System. The 7500 software uses the ROI calibration data to associate increases in fluorescence during a run with specific wells of the plate. The instrument uses a set of optic filters to distinguish the fluorescence emissions gathered during runs. You must generate a calibration image for each individual filter to account for minor differences in the optical path.

Note: The ROI calibration is a user-performed maintenance procedure.

Time Required 30 min

Materials Required



Applied Biosystems
Real-Time PCR
System Spectral
Calibration Kit



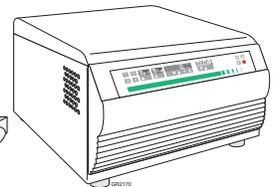
ROI Calibration
Plate



Powder-free
gloves



Safety
goggles



Centrifuge with
plate adapter

When to Perform the Calibration

Perform an ROI calibration:

- When installing the 7500/7500 Fast system. You must perform in sequence the ROI, background, optical, and dye calibrations and the instrument verification run.
- Every 6 months, or as often as necessary, depending on instrument use.
- After replacing the lamp.

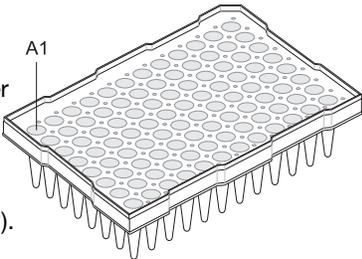
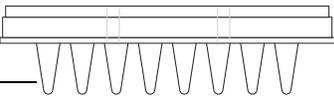
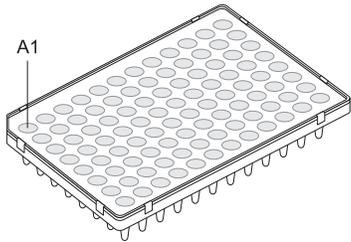
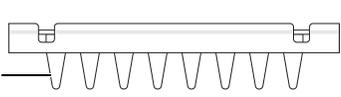
IMPORTANT! After every ROI calibration, you must perform a background calibration, optical calibration, dye calibration, and instrument verification.

Notes _____

Prepare the ROI Calibration Plate

Standard Plates Versus Fast Plates

Use the plate appropriate for your 7500/7500 Fast system.

Standard Plates – 7500 Systems	Fast Plates – 7500 Fast Systems
<p>Notched corner at top right</p>  <p>A1 at corner opposite from notched top-right corner (A12).</p>  <p>100-µL maximum reaction</p> <p>Vortex standard plates to ensure complete mixing, then centrifuge to ensure that all reagents are contained in the bottom of the well.</p> <p>CAUTION Optical 96-Well Fast Plates do not fit correctly into the standard block.</p>	<p>Notched corner at top left</p>  <p>A1 at notched top-left corner</p>  <p>30-µL maximum reaction volume</p> <p>Centrifuge Fast plates to ensure that all reagents are in the bottom of the well.</p> <p>Do not vortex Fast plates. Vortexing can cause loss of smaller reaction volumes.</p> <p>CAUTION Standard plates are not compatible with the 7500 Fast system and may be crushed by the 96-Well Fast Block.</p>

Prepare the Plate

IMPORTANT! Wear powder-free gloves when you handle the ROI calibration plate.

1. Obtain the ROI calibration plate from the spectral calibration kit in the freezer.
2. Allow the ROI calibration plate to warm to room temperature (approximately 5 min).

IMPORTANT! Do not remove an ROI calibration plate from its packaging until you are ready to run it. The fluorescent dye in the wells of the plate is photosensitive. Prolonged exposure to light can diminish the fluorescence from the plate.

Notes _____

- Remove the ROI calibration plate from its packaging. Leave the optical film on the plate.

IMPORTANT! Do not discard the packaging for the ROI calibration plate. The plate can be used up to three times if it is stored in its original packaging sleeve.



- (Standard plates only) Vortex the ROI calibration plate for 5 sec.

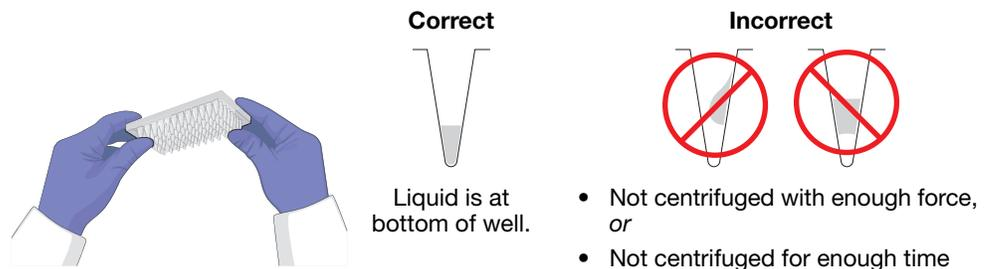
IMPORTANT! Do not vortex Fast plates.

(The remaining steps apply to both standard and Fast plates.)

- Centrifuge the plate for 2 min at less than 1500 rpm.

IMPORTANT! The ROI calibration plate must be well mixed and centrifuged.

- Verify that the liquid in each well of the ROI calibration plate is at the bottom of the well. If not, centrifuge the plate again at a higher rpm and for a longer period of time.



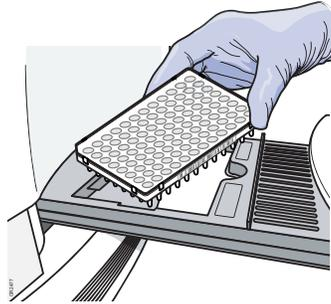
Notes _____

Load the Plate



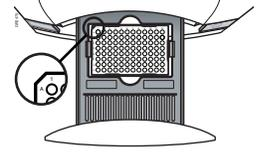
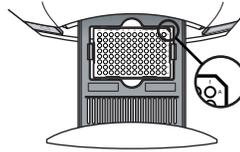
WARNING PHYSICAL INJURY HAZARD. During instrument operation, the sample block can be heated to 100 °C. Before performing the following procedure, be sure to wait until the sample block reaches room temperature.

1. Push the tray door to open it.
2. Load the plate into the plate holder in the instrument. Ensure that the plate is properly aligned in the holder.

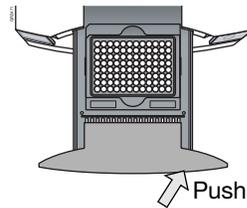
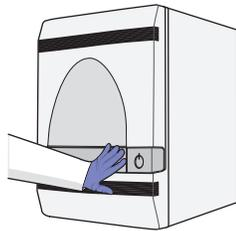


Load standard plates (7500 system) with the notched A12 position at the top-right of the tray.

Load Fast plates (7500 Fast system) with the notched A1 position at the top-left of the tray.



3. Close the tray door. Apply pressure to the right side of the tray door at an angle.



Notes

Perform the Calibration

Automated Versus Manual Calibrations

The 7500 software allows you to perform the ROI calibration automatically or manually.

ROI Calibration	Description	Use for...	See Page
Automatic	(Novice users) A wizard interface guides you through the calibration.	Regular calibrations	12
Manual	(Advanced users) An interface similar to previous versions of the software that allows you to perform the calibration manually.	<ul style="list-style-type: none"> Regular calibrations Custom dye calibration (see page 75) 	13

Perform an Automated ROI Calibration

Start the Calibration

1. In the 7500 software, select **Instrument ► Instrument Maintenance Manager**.
2. In the ROI screen of the Instrument Maintenance Manager, click **Start Calibration**.
3. Complete the calibration as instructed by the wizard.

The ROI Calibration dialog box displays three tabs:

- **Setup** – Displays instructions for setting up the ROI calibration. Clicking Next prompts opens the Run tab.
- **Run** – Clicking START RUN starts the calibration process and displays the processing messages. Clicking Next opens the Analysis tab.
- **Analysis** – Indicates the calibration status (Passed/Failed).

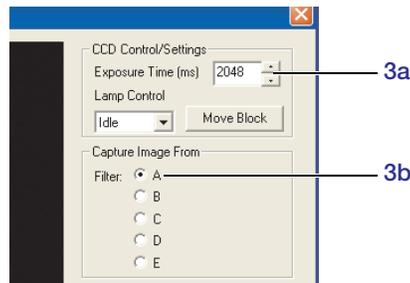
If you cannot obtain a passing calibration, see “[Troubleshoot the ROI Calibration](#)” on [page 17](#).

Notes _____

Perform a Manual ROI Calibration

Start the Calibration

1. In the 7500 software, select **Instrument ► Instrument Maintenance Manager**.
2. In the ROI tab of the Instrument Maintenance Manager, click **Start Manual Calibration**.
3. In the ROI Inspector dialog box:
 - a. In the Exposure Time field, enter **2048**.
 - b. Select **Filter A** (Filter 1).



4. Click **Snapshot** to generate an ROI image.
5. Determine if your ROI image is acceptable (the figures below show unsaturated and oversaturated images). Wells in an acceptable image:
 - Must be as bright as possible without oversaturating. (When you generate the ROI calibration, a warning is displayed if wells are oversaturated.)
 - Can contain some, but do not have to contain any, red pixels, which represent saturation.

Unsaturated	Normal Saturation	Over Saturated
Good	Acceptable	Not Acceptable

6. If your ROI image is acceptable, go to [step 7](#).

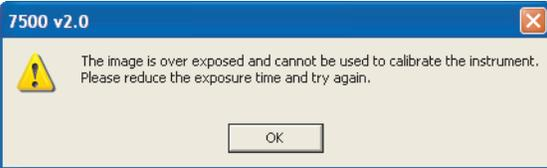
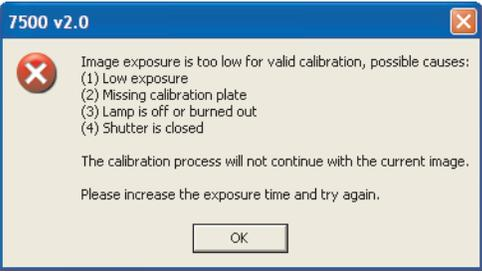
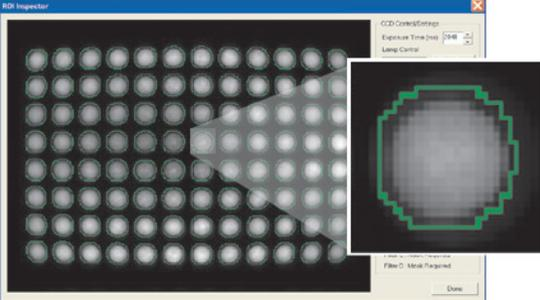
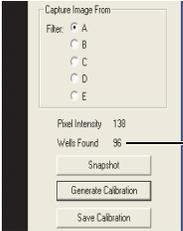
If your ROI image is oversaturated, decrease the Exposure Time by half, then click **Snapshot**. Repeat until you obtain an acceptable ROI image.

If you cannot obtain an acceptable image, see [“Troubleshoot the ROI Calibration” on page 17](#).

Notes

7. Click **Generate Calibration**.

The 7500 software takes a snapshot, then displays a message box or an ROI image:

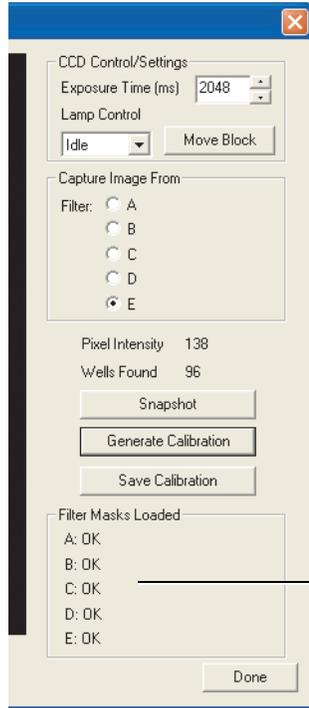
If the software displays:	Do the following:
 <p style="text-align: center;">or</p> 	<ol style="list-style-type: none"> Click OK. Decrease the Exposure Time by half (see step 3a on page 13). Repeat steps 3 and 5.
<p>A very faint ROI image, or you cannot generate a successful calibration.</p>	<p>See “Troubleshoot the ROI Calibration” on page 17.</p>
<p>An acceptable Calibration Image</p> <p>Note: Green circles appear around all wells indicating that the wells calibrated successfully.</p>  <p>Note: A green circle indicates that the region of interest for the well position is bright enough.</p>	<p>Verify that the:</p> <ul style="list-style-type: none"> Wells Found value is 96.  <ul style="list-style-type: none"> ROI Image displays a green circle around each well area.

8. Click **Save Calibration**.

The software saves the newly generated ROI calibration for Filter 1. “OK” appears next to Filter 1 in the Filter Masks Loaded section of the ROI Inspector.

Notes _____

- Repeat steps 3 through 8 for the remaining filters, resetting the Exposure Time to **2048** before performing the calibration for each filter. The ROI calibration is complete when Filter Masks Loaded for all the filters displays **OK**.



Calibration is complete when Filter Masks Loaded is **OK** for all Filters

- Click **Done**.

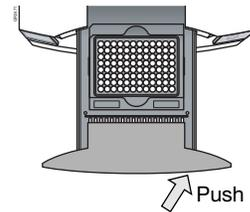
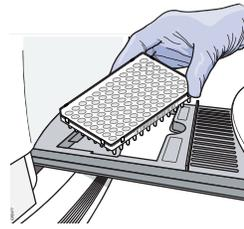
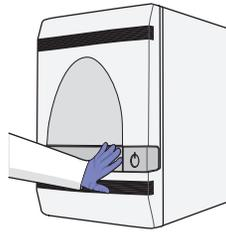
Notes

Unload the Plate



WARNING **PHYSICAL INJURY HAZARD.** During instrument operation, the sample block can be heated to 100 °C. Before performing the following procedure, be sure to wait until the sample block reaches room temperature.

1. Remove the calibration plate:
 - a. Push the tray door to open it.
 - b. Remove the calibration plate.
 - c. Push the tray door to close it.



2. Place the calibration plate inside its packaging sleeve. If you plan to perform background and optical calibrations:
 - Within the next 8 hr, keep the ROI calibration plate at room temperature. The optical calibration uses the ROI calibration plate.
 - On another day, return the packaged plate to the spectral calibration kit in the freezer.

IMPORTANT! Do not discard the calibration plate. If the plate is stored in its original packaging sleeve, you can use it up to three times after you open it.

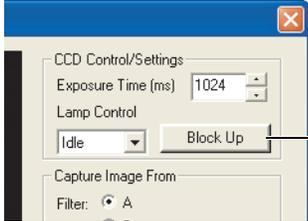
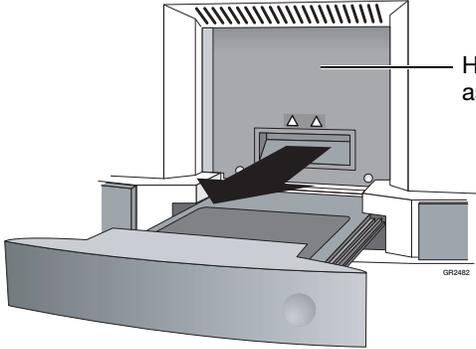


Continue with ““Perform the Background Calibration and Optical Calibration” on page 19.

IMPORTANT! After you perform an ROI calibration, you must also perform a background calibration (see page 20), an optical calibration (see page 25), dye calibrations (see page 32), and instrument verification (see page 46).

Notes _____

Troubleshoot the ROI Calibration

Problem/Symptom	Possible Cause	Action
ROI Calibration Failed	The sample block may be in its lowered position.	<ol style="list-style-type: none"> <li data-bbox="743 342 1382 396">1. If the CCD Control Settings in the ROI Inspector displays “Block Up,” click Block Up, to raise the block. <div data-bbox="776 411 1333 632">  </div>
ROI Image is Faint		<ol style="list-style-type: none"> <li data-bbox="743 646 1484 751">2. Check that the heated cover assembly is pulled all the way forward to ensure that the tray can be pushed in properly. If the 7500/7500 Fast system has a heated cover latch installed, check that the latch is in a locked position. <div data-bbox="776 766 1382 1115">  </div> <li data-bbox="743 1129 1484 1234">3. If the ROI calibration continues to fail, check the status of the halogen lamp within the 7500/7500 Fast instrument (see “Monitor the Lamp Status” on page 58), then replace the lamp if necessary (see “Replace the Halogen Lamp” on page 63).

Notes _____

Notes _____

3

Perform the Background Calibration and Optical Calibration

This chapter covers:

■ Perform the Background Calibration	20
■ Prepare the Background Calibration Plate	21
■ Perform the Background Calibration	23
■ Perform the Optical Calibration	25
■ Prepare the Calibration Plate	25
■ Perform the Optical Calibration	28
■ Troubleshoot the Background Calibration	30

Note: For more information about any of the topics discussed in this guide, access the Help from within Applied Biosystems 7500/7500 Fast Real-Time PCR System Software v2.0 by pressing **F1**, clicking  in the toolbar, or selecting **Help ▶ 7500 Software Help**.

Notes _____

Perform the Background Calibration

During a background calibration, the 7500/7500 Fast system:

- Performs reads of a background plate containing PCR buffer for 10 min at 60 °C.
- Averages the spectra recorded during the run and extracts the resulting spectral component to a calibration file.

The 7500 software then uses the calibration file during subsequent runs to remove the background fluorescence from the run data.

Note: The background calibration is a user-performed maintenance procedure.

Time Required 30 min

Materials Required



Applied Biosystems 7500/7500
Fast Real-Time PCR System
Spectral Calibration Kit

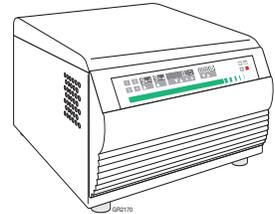
Background
Plate



Powder-free
gloves



Safety
goggles



Centrifuge with plate
adapter

When to Perform the Calibration

Perform a background calibration:

- When installing the 7500/7500 Fast system. You must perform in sequence the ROI, background, optical, and dye calibrations and the instrument verification run.
- Monthly, or as often as necessary, depending on instrument use.
- After replacing the lamp.

Background Fluorescence

Fluorescence data collected by the 7500/7500 Fast system includes a fluorescence signal inherent to the system, referred to as background fluorescence. Background fluorescence is a composite signal found in all spectral data. This signal consists of fluorescence from several sources, including:

- Background electronic signal
- Contaminants in the sample block
- The plastic consumable (plates and caps)

Guidelines for Calibration

- Make sure the centrifuge you use is clean. Before centrifuging, wipe down the bucket using a tissue.
- Handle the calibration plates with care to prevent contamination. Do not place plates on a lab bench, which may contaminate the plate. Always put calibration plates back into their original bags.

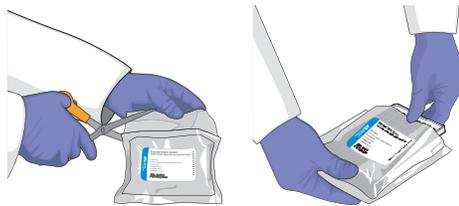
Notes _____

Prepare the Background Calibration Plate

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

1. Obtain the prepared background plate from the spectral calibration kit in the freezer.
2. Allow the background plate to warm to room temperature (at least 5 min).
3. Remove the background plate from its packaging.

IMPORTANT! Do not discard the packaging for the plate. The background plate can be used up to three times if it is stored in its original packaging sleeve.



4. (Standard plates only) Vortex the plate for 5 sec.

IMPORTANT! Do not vortex Fast plates.

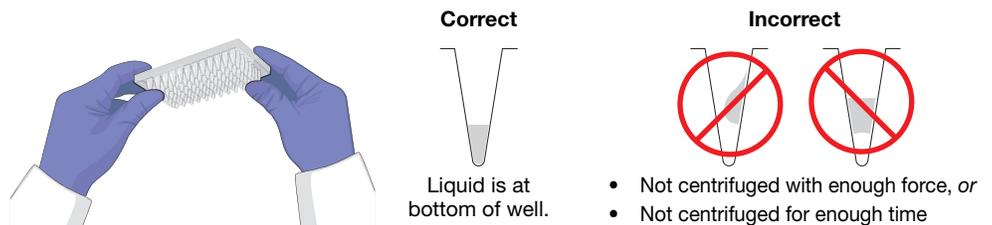
(The remaining steps apply to both standard and Fast plates.)

5. Centrifuge the plate for 2 min at less than 1500 rpm.

IMPORTANT! The plate must be well mixed and centrifuged.

6. Verify that the liquid in each well of the background plate is 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.



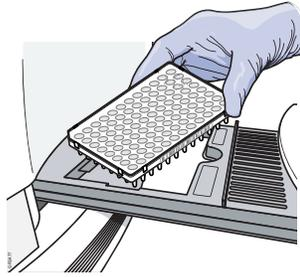
Notes

Load the Plate



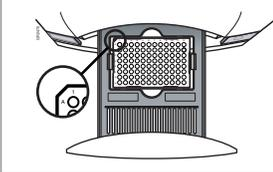
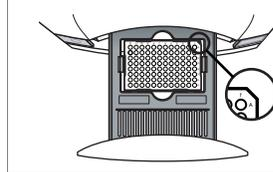
WARNING PHYSICAL INJURY HAZARD. During instrument operation, the sample block can be heated to 100 °C. Before performing the following procedure, be sure to wait until the sample block reaches room temperature.

1. Push the tray door to open it.
2. Load the plate into the plate holder in the instrument. Ensure that the plate is properly aligned in the holder.

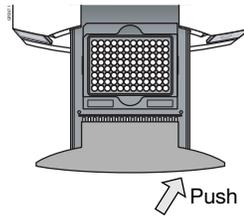
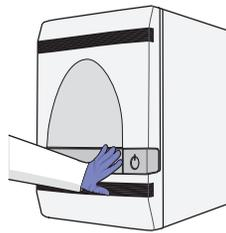


Load standard plates (7500 system) with the notched A12 position at the top-right of the tray.

Load Fast plates (7500 Fast system) with the notched A1 position at the top-left of the tray.



3. Close the tray door. Apply pressure to the right side of the tray door at an angle.



Note: If you cannot open the tray, the sample block may be in its raised position, locking the tray door position. To lower the block, select **Instrument ▶ Calibrate**, then exit the ROI Inspector.

Notes _____

Perform the Background Calibration

Perform the Calibration

1. In the 7500 software, select **Instrument ▶ Instrument Maintenance Manager**.
2. In the Instrument Maintenance Manager, select the **Background** tab.
3. In the Background tab, click **Start Calibration**.
4. Complete the calibration as instructed by the wizard.

The Background Calibration dialog box displays four tabs:

- **Overview** – Displays information describing the calibration.
- **Setup** – Displays instructions for setting up the background calibration. Clicking Next prompts opens the Run tab.
- **Run** – Clicking START RUN starts the calibration process and displays the processing messages. Clicking Next opens the Analysis tab.
- **Analysis** – Indicates the calibration status (Passed/Failed).

If you cannot obtain a passing calibration, see [“Troubleshoot the Background Calibration”](#) on page 30.

Note: Before starting the calibration, the instrument may pause (up to 10 min) to allow the heated cover to reach temperature.

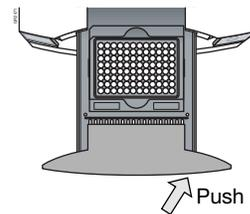
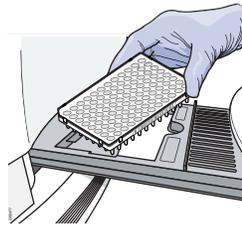
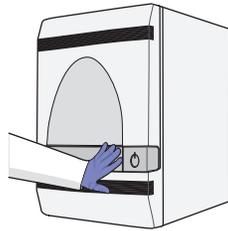
Notes _____

Unload the Plate



WARNING PHYSICAL INJURY HAZARD. During instrument operation, the sample block can be heated to 100 °C. Before performing the following procedure, be sure to wait until the sample block reaches room temperature.

1. Remove the calibration plate:
 - a. Push the tray door to open it.
 - b. Remove the calibration plate.
 - c. Push the tray door to move it into the instrument.



2. Place the calibration plate inside its packaging sleeve, then return the packaged plate to the spectral calibration kit in the freezer.

IMPORTANT! Do not discard the calibration plate. If the plate is stored in its original packaging sleeve, you can use the plate up to three times after you open it.



If necessary, continue with “[Perform the Optical Calibration](#)” on page 25.

You must perform an optical calibration (see [page 25](#)), dye calibrations (see [page 32](#)), and instrument verification (see [page 46](#)) if you are performing the background calibration:

- As part of your semiannual maintenance
- After replacing or moving any parts of the optics

For more information, see “[Recommended Maintenance Schedule](#)” on page 5.

Notes _____

Perform the Optical Calibration

The optical calibration generates data that allows the 7500 software to compensate for the physical effects of the fifth filter in the 7500/7500 Fast system.

Note: The optical calibration is a user-performed maintenance procedure.

Time Required 10 min

Materials Required ROI calibration plate

When to perform the Calibration Perform an optical calibration:

- When installing the 7500/7500 Fast system. You must perform in sequence the ROI, background, optical, and dye calibrations and the instrument verification run.
- Every 6 months, or as often as necessary, depending on instrument use.

Prepare the Calibration Plate

Prepare the Plate If you kept your ROI calibration plate at room temperature after performing an ROI calibration (see [Chapter 2](#)), skip to [step 5 on page 26](#) to spin down any condensation that may have formed when the plate was at room temperature. If the ROI calibration plate is in the freezer, go to [step 1](#).

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

1. Obtain the ROI calibration plate from the spectral calibration kit in the freezer.
2. Allow the ROI calibration plate to warm to room temperature (at least 5 min).
3. Remove the ROI calibration plate from its packaging.

IMPORTANT! Do not discard the packaging for the plate. The ROI calibration plate can be used up to three times if it is stored in its original packaging sleeve.



Notes _____

4. (Standard plates only) Vortex the plate for 5 sec.

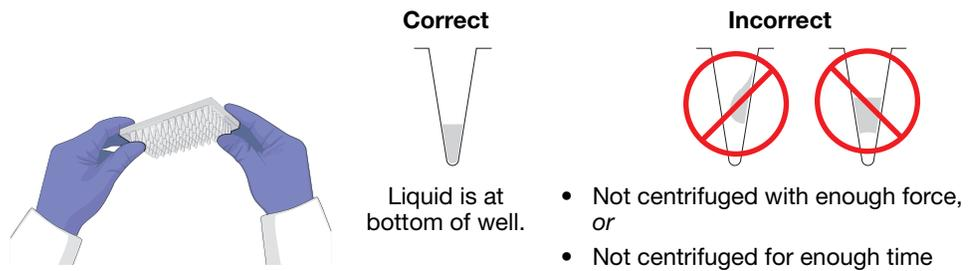
IMPORTANT! Do not vortex Fast plates.

(The remaining steps apply to both standard and Fast plates.)

5. Centrifuge the plate for 2 min at less than 1500 rpm.

IMPORTANT! The ROI calibration plate must be well mixed and centrifuged.

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



Notes _____

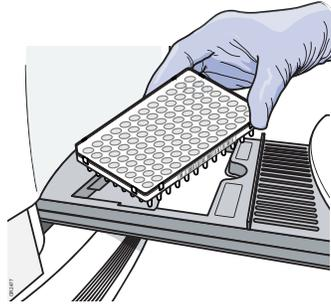
Load the Plate



WARNING

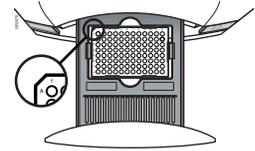
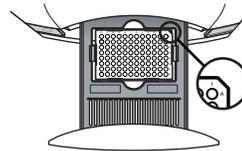
PHYSICAL INJURY HAZARD. During instrument operation, the sample block can be heated to 100 °C. Before performing the following procedure, be sure to wait until the sample block reaches room temperature.

1. Push the tray door to open it.
2. Load the plate into the plate holder in the instrument. Ensure that the plate is properly aligned in the holder.

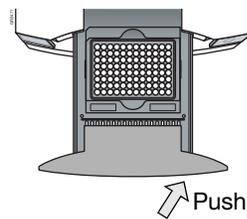
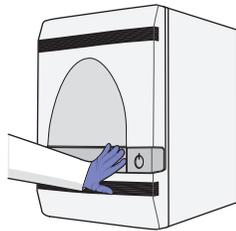


Load standard plates (7500 system) with the notched A12 position at the top-right of the tray.

Load Fast plates (7500 Fast system) with the notched A1 position at the top-left of the tray.



3. Close the tray door. Apply pressure to the right side of the tray door at an angle.



Notes _____

Perform the Optical Calibration

Perform the Calibration

1. In the 7500 software, select **Instrument ▶ Instrument Maintenance Manager**.
2. In the Instrument Maintenance Manager, select the **Optical** tab.
3. In the Optical screen, click **Start Calibration**.
4. Complete the calibration as instructed by the wizard.

The Optical Calibration dialog box displays four tabs:

- **Overview** – Displays information describing the calibration.
- **Setup** – Displays instructions for setting up the optical calibration. Clicking Next prompts opens the Run tab.
- **Run** – Clicking START RUN starts the calibration process and displays the processing messages. Clicking Next opens the Analysis tab.
- **Analysis** – Indicates the calibration status (Passed/Failed).

If you cannot obtain a passing calibration, see [“Troubleshoot the Background Calibration”](#) on page 30.

Note: Before starting the calibration, the instrument may pause (up to 10 min) to allow the heated cover to reach temperature.

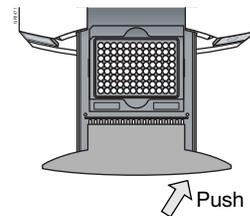
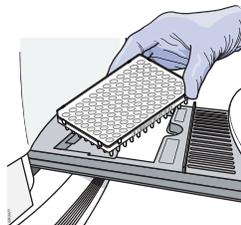
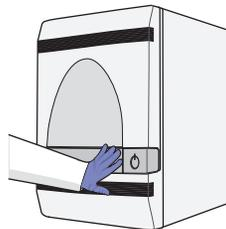
Unload the Plate



WARNING

PHYSICAL INJURY HAZARD. During instrument operation, the sample block can be heated to 100 °C. Before performing the following procedure, be sure to wait until the sample block reaches room temperature.

1. Remove the calibration plate:
 - a. Push the tray door to open it.
 - b. Remove the calibration plate.
 - c. Push the tray door to move it into the instrument.



Notes _____

2. Place the calibration plate inside its packaging sleeve. Return the packaged plate to the spectral calibration kit in the freezer.

IMPORTANT! Do not discard the calibration plate. If the plate is stored in its original packaging sleeve, you can use it up to three times after you open it.



Continue with “Perform the Dye Calibration” on page 31.

Notes _____

Troubleshoot the Background Calibration

Problem/Symptom	Possible Cause	Action
Background Calibration Failed	One or more wells of the background plate produced spectra that exceed the maximum limit for the instrument.	<ol style="list-style-type: none"> 1. Repeat the calibration using the same background plate. 2. If the calibration fails again, repeat the calibration using a different background plate. 3. If the calibration fails again, determine the source of the contamination, as explained in “How to Identify Contamination” below.

How to Identify Contamination

Signals that exceed the limit of normal background fluorescence may indicate fluorescent contaminants on the calibration plate or the sample block. Common contaminants include ink residue from permanent pens, powder from disposable gloves, and dust.

To determine the source and location of the contamination:

1. While viewing the raw spectra, locate the contaminated well position(s) by selecting successively smaller regions of the plate layout.
2. Rotate the background plate 180°, then perform the background calibration again.
3. Repeat [step 1](#) to locate the contamination. If the well position(s) of the contamination in [steps 1](#) and [3](#) are:
 - **Identical** – The sample block is contaminated. Decontaminate the sample block as explained in “[Decontaminate the Sample Block](#)” on [page 59](#).
 - **Reversed** – The background plate is contaminated. Discard the background plate and perform the background run using a new background plate.

If the calibration fails after you replace the background plate and decontaminate the sample block:

1. Load a plate covered by a piece of black paper into the 7500/7500 Fast instrument.
2. Perform the background run as explained in this chapter.
3. After the run is complete, select all wells of the plate layout.
4. View the Spectral plot for the peak(s). If a peak is:
 - **Visible** – The optics of your 7500/7500 Fast instrument may be contaminated. Contact Applied Biosystems as explained in “[How to Obtain Support](#)” on [page x](#).
 - **Absent** – The sample block is contaminated. Decontaminate the sample block as explained in “[Decontaminate the Sample Block](#)” on [page 59](#).

Notes _____

4

Perform the Dye Calibration

This chapter covers:

- Overview 32
- Prepare the Dye Calibration Plates 35
- Perform the Dye Calibration 36
- Troubleshoot the Dye Calibration 43

Note: For more information about any of the topics discussed in this guide, access the Help from within Applied Biosystems 7500/7500 Fast Real-Time PCR System Software v2.0 by pressing **F1**, clicking  in the toolbar, or selecting **Help ▶ 7500 Software Help**.

Notes

Overview

During a dye calibration, the Applied Biosystems 7500/7500 Fast Real-Time PCR System:

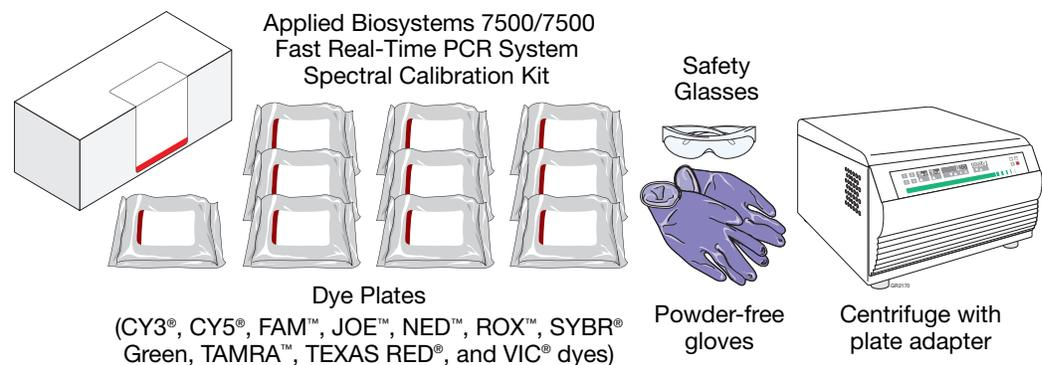
- Collects spectral data from a series of dye standards.
- Stores the spectral information for the dye standards in a pure spectra calibration file.

The software uses the pure spectra data during experiment runs to characterize and distinguish the individual contribution of each dye in the total fluorescence collected by the instrument. After each run, the 7500 software receives data in the form of a raw spectra signal for each reading. It determines the contribution of each fluorescent dye used in the sample by comparing the raw spectra to the pure spectra calibration data. When you save an experiment after analysis, the software stores the pure spectra with the collected fluorescence data for that experiment.

Note: The dye calibration is a user-performed maintenance procedure.

Time Required 1 hr

Materials Required



Note: If you store Applied Biosystems 7500/7500 Fast Real-Time PCR System dye plates in their original packaging in the freezer, you can use them to calibrate a 7500/7500 Fast instrument up to 3 times for 6 months after opening them.

When to Perform Dye Calibration

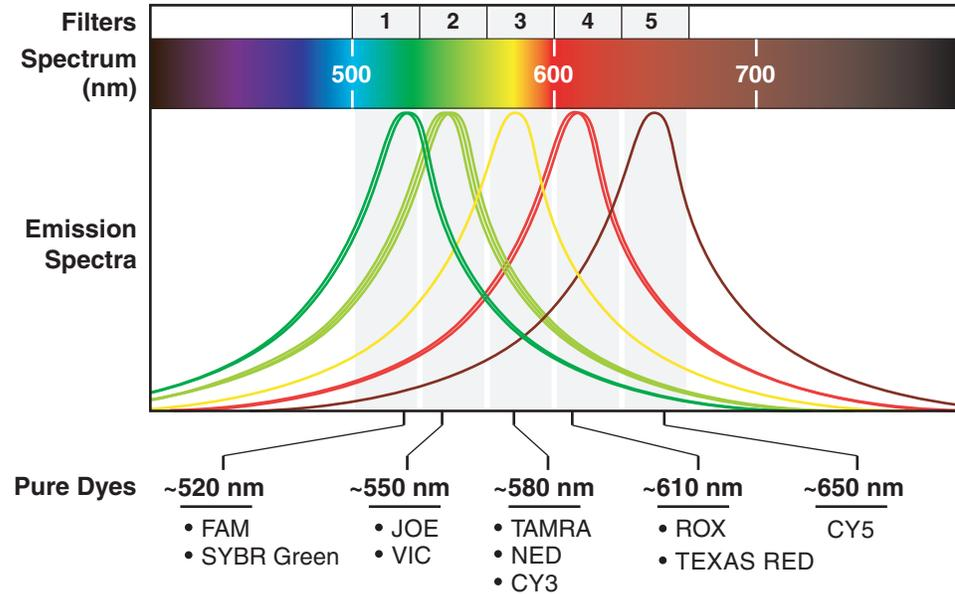
Perform a dye calibration:

- When installing the 7500/7500 Fast system. You must perform in sequence the ROI, background, optical, and dye calibrations and the instrument verification run.
- Every 6 months, or as often as necessary, depending on instrument use.

IMPORTANT! You must perform a background run before every series of dye calibrations. Because the age and use of instrument components can affect pure spectra readings, Applied Biosystems recommends performing a dye calibration at least every 6 months.

Notes _____

Dye Sets The Applied Biosystems 7500/7500 Fast Real-Time PCR Systems use the following dye sets for calibration: CY3[®] dye, CY5[®] dye, FAM[™] dye, JOE[™] dye, NED[™] dye, ROX[™] dye, SYBR[®] Green dye, TAMRA[™] dye, TEXAS RED[®] dye, and VIC[®] dye. The following figure shows the emission spectrum for each dye, and the filters and wavelengths at which each dye is read.



Custom Dye The 7500/7500 Fast system can be used to run assays designed with custom dyes (dyes not supplied by Applied Biosystems). However, before using custom dyes with the 7500/7500 Fast instrument, you must create and run a custom calibration plate. The 7500 software uses the custom calibration plate to create a spectral standard to distinguish the custom dye in the fluorescence data collected during the run. See [Appendix B](#) for information on custom dye calibrations.

IMPORTANT! To use a custom dye on your 7500/7500 Fast system, it must fluoresce within the 520 to 650 nm spectral range measured by the 7500/7500 Fast instrument.

Notes _____

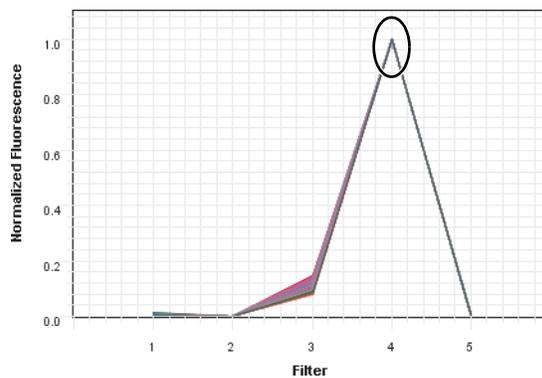
About the Analysis

The product of a dye calibration is a collection of spectral profiles that represent the fluorescence signature of each dye standard. Each profile consists of a set of spectra that correspond to the fluorescence collected from the wells of the spectral calibration plate. The 7500 software plots the resulting data for each spectral profile in a graph of fluorescence versus filter.

When the 7500 software extracts the calibration data from a dye run, it evaluates the fluorescence signal generated by each well in terms of the collective spectra for the entire calibration plate. Dye spectra are generally acceptable if they peak within the same filter but diverge slightly at other wavelengths (see below).

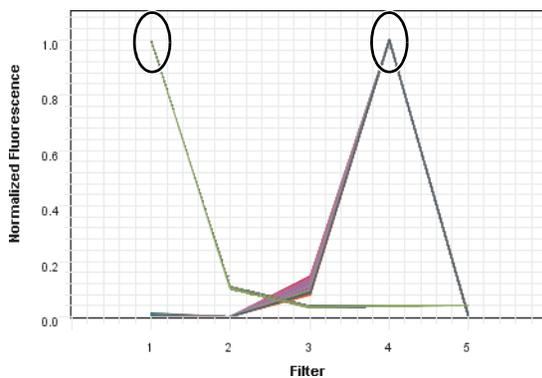
The 7500 software can compensate for some differences in a spectral profile by replacing (auto-repairing) the spectra of unacceptable wells with the spectra of other wells on the reaction plate. However, the software allows only a few replacements and may reject the calibration if the spectra between neighboring wells vary significantly.

Note: Because the wells in a calibration plate contain dyes at identical concentrations, the resulting signals for the wells containing each dye should be similar. Among wells containing the same dye, variations in spectral position and peak position are caused by minor differences in the optical properties and excitation energy between the individual wells.



Acceptable Spectra

Spectra peak at the same wavelength and do not diverge significantly



Unacceptable Spectra

Spectra peak at the different wavelengths

Notes

Prepare the Dye Calibration Plates

IMPORTANT! Before performing a dye calibration, you must perform an ROI calibration (see [page 8](#)), a background calibration (see [page 20](#)), and an optical calibration (see [page 25](#)).

Prepare the Plates

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

1. Obtain the spectral calibration kit from the freezer, then remove all of the dye plates.
 2. Return the spectral calibration kit to the freezer.
 3. Allow the dye plates to warm to room temperature (approximately 5 min).
-

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

Note: If you store Applied Biosystems 7500/7500 Fast Real-Time PCR System dye plates in their original packaging in the freezer, you can use them to calibrate a 7500/7500 Fast instrument up to 3 times for 6 months after opening them.

Continue with “[Perform the Dye Calibration](#)” on [page 36](#).

Notes _____

Perform the Dye Calibration

Perform the Calibration

1. In the 7500 software, select **Instrument ► Instrument Maintenance Manager**.
2. In the Instrument Maintenance Manager, select the **Dye** tab.
3. In the Dye screen, select **System Dye Calibration**.
4. Click **Start Calibration**.
5. Complete the calibration for each plate as instructed by the wizard.

IMPORTANT! The wizard guides you through the calibration of each dye separately. You must set up, run, and analyze each dye plate independently.

The Dye Calibration dialog box displays four tabs:

- **Overview** – Displays information describing the calibration.
When the software prompts you to obtain the required materials, select the dyes that you want to calibrate.
- **Setup** – Displays instructions for setting up the dye calibration. Clicking Next prompts opens the Run tab.
When the software prompts you to load each dye plate, prepare and load the plates as described in [“Load a Dye Plate” on page 37](#).
- **Run** – Clicking START RUN starts the calibration process and displays the processing messages. Clicking Next opens the Analysis tab.
- **Analysis** – Indicates the calibration status (Passed/Failed).

When the software prompts you to analyze the spectra collected from each dye plate, verify the status of the calibration:

- **Passed** – The 7500/7500 Fast instrument passed the calibration. Go to [“Analyze the Calibration Data” on page 39](#).
- **Failed** – The 7500/7500 Fast instrument failed the calibration. Troubleshoot the error as described in [“Troubleshoot the Dye Calibration” on page 43](#).

Note: Before starting the calibration, the instrument may pause (up to 10 min) to allow the heated cover to reach temperature.

Load a Dye Plate

Note: Because the wizard guides you through the calibration of each dye separately, perform the following procedure for each dye that you calibrate.

1. Remove the dye plate that is specified by the software from its packaging.

IMPORTANT! Do not discard the packaging for the plate. The plate can be used up to three times if it is stored in its original packaging sleeve.



2. (Standard plates only) Vortex the plate for 5 sec.

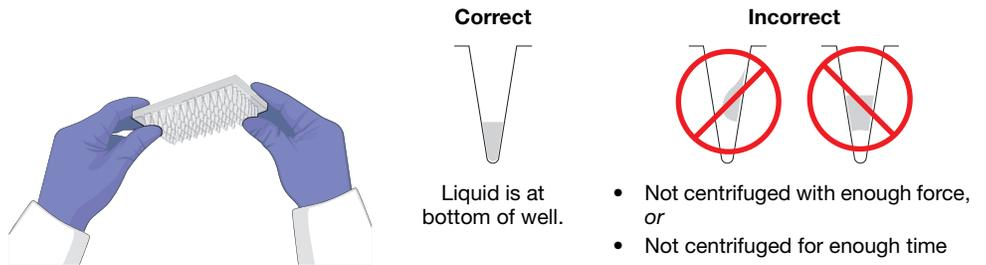
IMPORTANT! Do not vortex Fast plates.

(The remaining steps apply to both standard and Fast plates.)

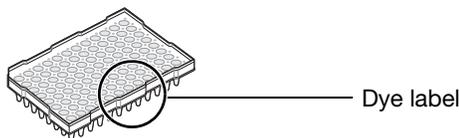
3. Centrifuge the plate for 2 min at less than 1500 rpm.

IMPORTANT! The plate must be well mixed and centrifuged.

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



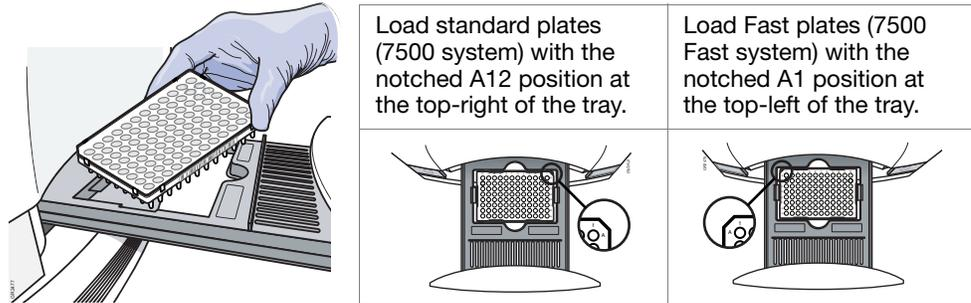
5. Verify that the dye plate that you are about to load matches the dye selected in the 7500 software.



6. Push the tray door to open it.

Notes

7. Load the plate into the plate holder in the instrument. Ensure that the plate is properly aligned in the holder.



8. Close the tray door. Apply pressure to the right side of the tray door at an angle.

Note: If you cannot open the tray, the sample block may be in its raised position, locking the tray position. To lower the block, select **Instrument ▶ Calibrate**, then exit the ROI Inspector.

Notes _____

Analyze the Calibration Data

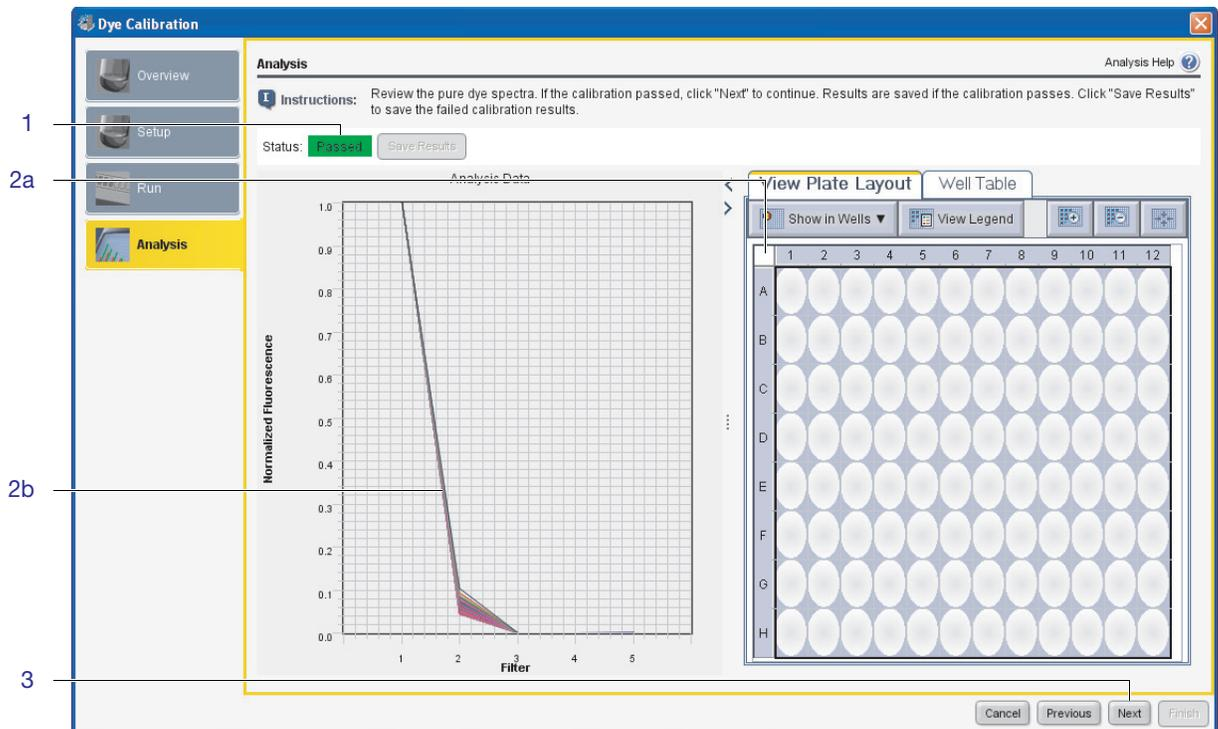
Note: Because the wizard guides you through the calibration of each dye separately, perform the following procedure for each dye that you calibrate.

1. Verify the status of the calibration:
 - **Passed** – The 7500/7500 Fast instrument passed the calibration. Go to [step 2](#).
 - **Failed** – The 7500/7500 Fast instrument failed the calibration. Troubleshoot the error as described in [“Troubleshoot the Dye Calibration”](#) on page 43.
2. Verify the grouping of the dye spectra:
 - a. In the plate layout, select the wells of the plate.
 - b. Inspect the raw data. For each spectrum, verify that the peak is:
 - Within the detectable range for the 7500/7500 Fast instrument.
 - Free of irregular spectral peaks.
 - Present in the correct channel for the dye (see [Table 1](#) on page 41).

If a spectrum does not match the criteria above, troubleshoot the problem as described in [“Troubleshoot the Dye Calibration”](#) on page 43.

Note: Among wells containing the same dye, variations in spectral position and peak position are caused by minor differences in the optical properties and excitation energy between the individual wells.

3. If all spectra are acceptable, then click **Next**.

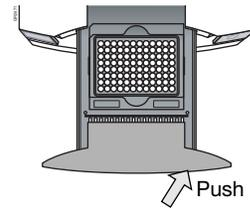
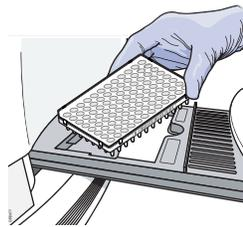
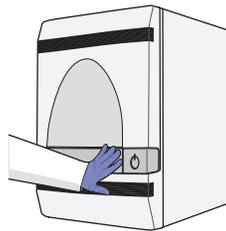


Notes

4. Remove the calibration plate:

WARNING PHYSICAL INJURY HAZARD. During instrument operation, the sample block can be heated to 100 °C. Before performing the following procedure, be sure to wait until the sample block reaches room temperature.

- a. Push the tray door to open it.
- b. Remove the calibration plate.
- c. Push the tray door to close it.



- d. Place the calibration plate inside its packaging sleeve. Return the packaged plate to the spectral calibration kit in the freezer.

Note: If you store Applied Biosystems 7500/7500 Fast Real-Time PCR System dye plates in their original packaging in the freezer, you can use them to calibrate 7500/7500 Fast instruments up to 3 times for 6 months after opening them.



5. After you remove the dye plate as instructed, click **Finish**.
6. Prepare and run the next plate as explained in “[Prepare the Plates](#)” on page 35.

Notes _____

Table 1 7500/7500 Fast system dye spectra

Filter	Peak (nm)	Dye/Spectra		
1	~520	<p>FAM dye</p>	<p>SYBR Green dye</p>	
2	~550	<p>JOE dye</p>	<p>VIC dye</p>	
3	~580	<p>CY3 dye</p>	<p>NED dye</p>	<p>TAMRA dye</p>

Notes

Table 1 7500/7500 Fast system dye spectra

Filter	Peak (nm)	Dye/Spectra	
4	~610	<p>ROX dye</p>	<p>TEXAS RED dye</p>
5	~670	<p>CY5 dye</p>	

Notes

Troubleshoot the Dye Calibration

Problem/Symptom	Possible Cause	Action
One or more raw spectra are at or below the detectable threshold for the calibration.	<ul style="list-style-type: none"> The spectral calibration plate was centrifuged insufficiently. The spectral calibration plate contains old or insufficient reagents. If you are running a custom spectral calibration plate, the dye may not be present at a sufficient concentration. 	<ol style="list-style-type: none"> Unload the 7500/7500 Fast instrument and view the wells of the spectral calibration plate. If the liquid in the wells is not: <ul style="list-style-type: none"> At the bottom of the wells, centrifuge the plate for a longer time, then repeat the calibration. Equivalent in volume, the plate is not sealed and the reagents have evaporated. Discard it and run another. If the spectral calibration plate appears to be normal, discard the plate and run another. If the problem persists, contact Applied Biosystems as explained in “How to Obtain Support” on page x. <p>Note: If you are running a custom spectral calibration plate, create another plate but increase the concentration of the dye that produced insufficient signal.</p>
One or more raw spectra exceed the maximum limit for the 7500/7500 Fast instrument.	<ul style="list-style-type: none"> Fluorescent contaminants are on the sample block(s) or spectral calibration plate. If you are running a custom spectral calibration plate, the dye may be too concentrated. 	<p>Verify that contaminants are not present by performing a background calibration as explained in Chapter 3, “Perform the Background Calibration and Optical Calibration.” If the background calibration does not show sample block contamination, the spectral calibration plate may be contaminated.</p> <p>Note: If you are running a custom spectral calibration plate, create another plate but decrease the concentration of the dye that exceeds the detectable limit.</p>
The spectra contain peaks in more than one filter.	Fluorescent contaminants are on the sample block(s) or spectral calibration plate.	

Notes _____

Notes _____

5

Verify the Instrument Performance

This chapter covers:

■ Overview.....	46
■ Set Up the Experiment.....	48
■ Run the Experiment.....	49
■ Analyze the Experiment.....	51
■ Troubleshoot the RNase P Experiment.....	54

Note: For more information about any of the topics discussed in this guide, access the Help from within Applied Biosystems 7500/7500 Fast Real-Time PCR System Software v2.0 by pressing **F1**, clicking  in the toolbar, or selecting **Help ▶ 7500 Software Help**.

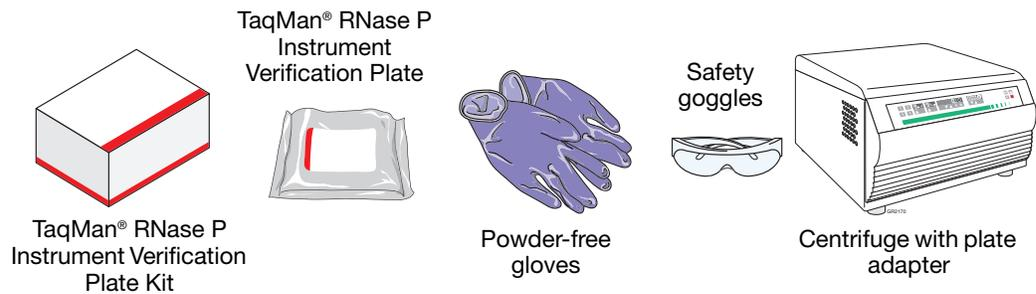
Notes

Overview

Perform the TaqMan® RNase P Instrument Verification Plate run to verify the performance of an Applied Biosystems 7500/7500 Fast Real-Time PCR System.

Time Required 1 hr

Materials Required



When to Perform the RNase P Experiment

Applied Biosystems recommends performing an RNase P experiment:

- When installing the 7500/7500 Fast system. You must perform in sequence the ROI, background, optical, and dye calibrations and the instrument verification run.
- After moving the instrument to another location.
- As needed to verify the function of the 7500/7500 Fast instrument.

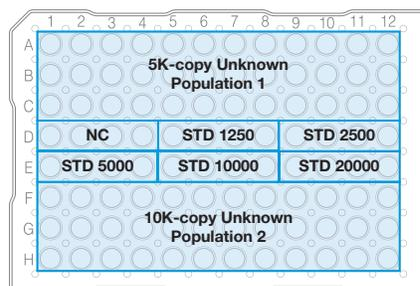
About the RNase P Plate

The RNase P plate is preloaded with the reagents necessary for the detection and quantitation of genomic copies of the human RNase P gene (a single-copy gene encoding the RNase moiety of the RNase P enzyme).

Each well contains:

- 1X TaqMan® Fast Universal PCR Master Mix, No AmpErase® UNG
- RNase P primers
- FAM™ dye-labeled probe
- Known concentration of human genomic DNA template

The figure below illustrates the arrangement of the standard and unknown populations on the RNase P plate. The RNase P plate contains five replicate groups of standards (1250, 2500, 5000, 10,000, and 20,000 copies), two unknown populations (5000 and 10,000 copies), and negative control wells (NC).



Notes _____

After the run, the 7500 software:

1. Generates a standard curve from the averaged threshold cycle (C_T) values of the replicate groups of standards.
2. Calculates the concentration of the two unknown populations using the standard curve.
3. Calculates the following to assess the 7500/7500 Fast instrument performance:

$$[(CopyUnk_2) - 3(\sigma_{CopyUnk_2})] > [(CopyUnk_1) + 3(\sigma_{CopyUnk_1})]$$

where:

- $CopyUnk_1$ = Average copy number of unknown #1 (5,000-copy population)
- $\sigma_{CopyUnk_1}$ = Standard deviation of unknown #1 (5,000-copy population)
- $CopyUnk_2$ = Average copy number of unknown #2 (10,000-copy population)
- $\sigma_{CopyUnk_2}$ = Standard deviation of unknown #2 (10,000-copy population)

Installation Specification

The 7500/7500 Fast instrument passes the installation specification if the inequality holds and the instrument successfully distinguishes between 5,000 and 10,000 copies with a statistical confidence level of 99.7%.

To meet the installation specification, you can omit a limited number of outlier wells from the 5,000- and 10,000-copy unknown populations.

Instrument	Maximum Number of Outlier Wells That Can Be Removed			
	Unknown Population		Standard (STD)	Negative Controls (NC)
	5,000-copy	10,000-copy		
7500 System	6	6	0	0
7500 Fast System				

Notes

Set Up the Experiment

Prepare the TaqMan® RNase P Fast Instrument Verification Plate for the run.

Prepare the RNase P Plate

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

1. Obtain the TaqMan® RNase P Fast Instrument Verification Plate from the freezer, then allow the reaction plate to warm to room temperature (for approximately 5 min).
2. Remove the RNase P plate from its packaging.



3. (Standard plates only) Vortex the plate for 5 sec.

IMPORTANT! Do not vortex Fast plates.

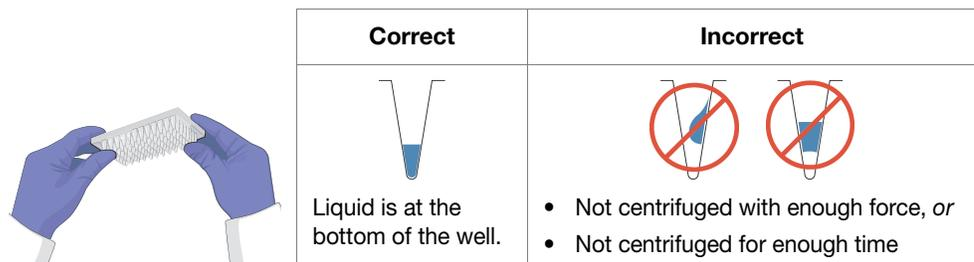
(The remaining steps apply to both standard and Fast plates.)

4. Centrifuge the reaction plate for 2 min at less than 1500 rpm.

IMPORTANT! The reaction plate must be well mixed and centrifuged.

5. Verify that the liquid is at the bottom of each well of the reaction plate. If not, centrifuge the reaction plate again at a greater rpm and for a longer time.

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



Notes _____

Run the Experiment

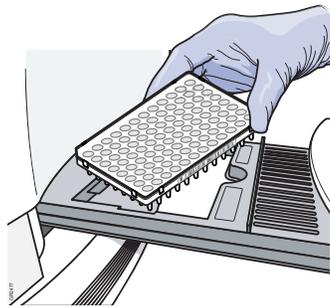
After preparing the TaqMan® RNase P Fast Instrument Verification Plate, load the plate into the 7500/7500 Fast instrument and start the run.

Load the Plate



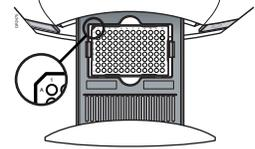
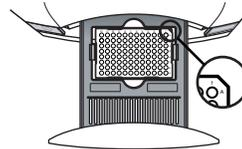
WARNING PHYSICAL INJURY HAZARD. During instrument operation, the sample block can be heated to 100 °C. Before performing the following procedure, be sure to wait until the sample block reaches room temperature.

1. Push the tray door to open it.
2. Load the plate into the plate holder in the instrument. Ensure that the plate is properly aligned in the holder.

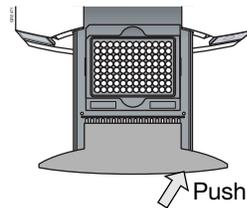
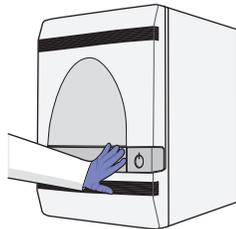


Load standard plates (7500 system) with the notched A12 position at the top-right of the tray.

Load Fast plates (7500 Fast system) with the notched A1 position at the top-left of the tray.



3. Close the tray door. Apply pressure to the right side of the tray door at an angle.



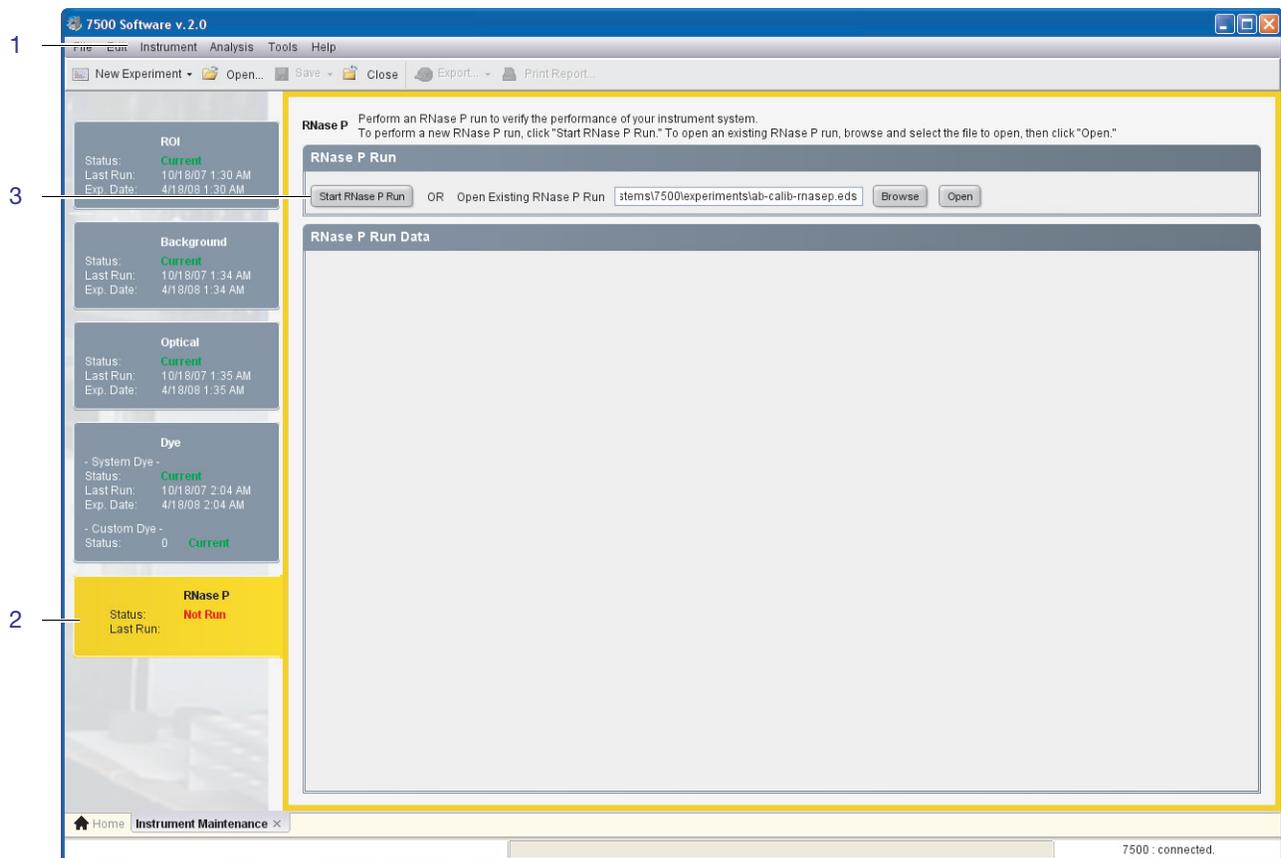
Notes _____

- Start the Run**
1. In the 7500 software, select **Instrument ► Instrument Maintenance Manager**.
 2. In the Instrument Maintenance Manager, select the **RNase P** tab.
 3. In the RNase P screen, click **Start RNase P Run**.
 4. Complete the calibration as instructed by the wizard.

The RNase P dialog box displays four tabs:

- **Overview** – Displays information describing the calibration.
- **Setup** – Displays instructions for setting up the RNase P run. Clicking Next prompts opens the Run tab.
- **Run** – Clicking START RUN starts the run process and displays the processing messages. Clicking Next opens the Analysis tab.
- **Analysis** – Indicates the run status (Passed/Failed).

Note: Before starting the run, the instrument may pause (up to 10 min) to allow the heated cover to reach the correct temperature.



Notes _____

Analyze the Experiment

Review the data to verify the results of the experiment.

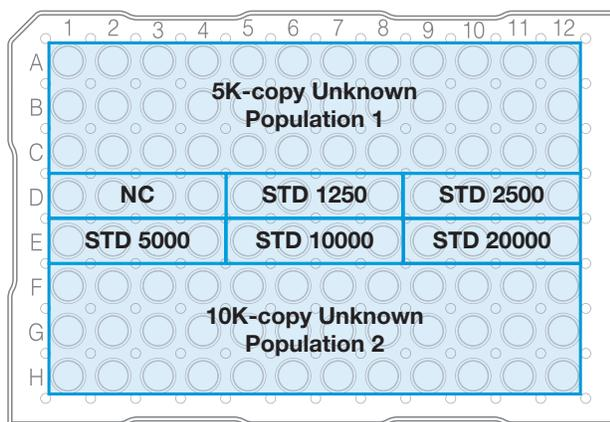
Verify the Results of the Analysis

Note: After the 7500 software completes the RNase P run, it automatically analyzes the run and displays the results in the Analysis screen.

1. In the Analysis screen of the RNase P Run wizard, verify the status of the run:
 - **Passed** – The 7500/7500 Fast instrument passed the RNase P run. Go to [step 5 on page 53](#).
 - **Failed** – The 7500/7500 Fast instrument failed the RNase P run. Go to [step 2](#) to review the data for outliers.

If the run fails, the automated analysis may have included outliers that caused the initial analysis to fail. Experimental error may cause some wells to be amplified insufficiently or not at all. These wells typically produce C_T values that differ significantly from the average for the associated replicate wells. If included in the calculations, these outlying data (outliers) can result in erroneous measurements.

2. In the Amplification Plot, select **Ct vs. Well** from the Plot Type drop-down list.
3. Verify the uniformity of each replicate population on the reaction plate (controls, standards, and unknowns) by comparing the groupings of C_T values:
 - a. In the plate layout, select the wells containing the 10,000-copy unknown population (wells rows F, G, and H).



- b. In the plot, verify that the C_T s of the replicate population are equivalent.

Note: The numbers on the X-axis of the plot correspond to the wells of the reaction plate. Beginning with well A1, the wells are numbered from left-to-right, and top-to-bottom.

Notes

- c. If an outlier is in the selected population, select the corresponding well in the plate layout, then click **Omit** to remove the well from the analysis.

Instrument	Maximum Number of Outlier Wells That Can Be Removed			
	Unknown Population		Standard (STD)	Negative Control (NC)
	5,000-copy	10,000-copy		
7500 System	6	6	0	0
7500 Fast System				

IMPORTANT! If the number of outliers exceeds the limit in the table above, order another RNase P plate and repeat the experiment.

- d. Repeat [steps 3a](#) through [3c](#) for each replicate population (unknowns, standards, and negative controls) on the reaction plate.
4. Click **Reanalyze** to analyze the run without the outliers.

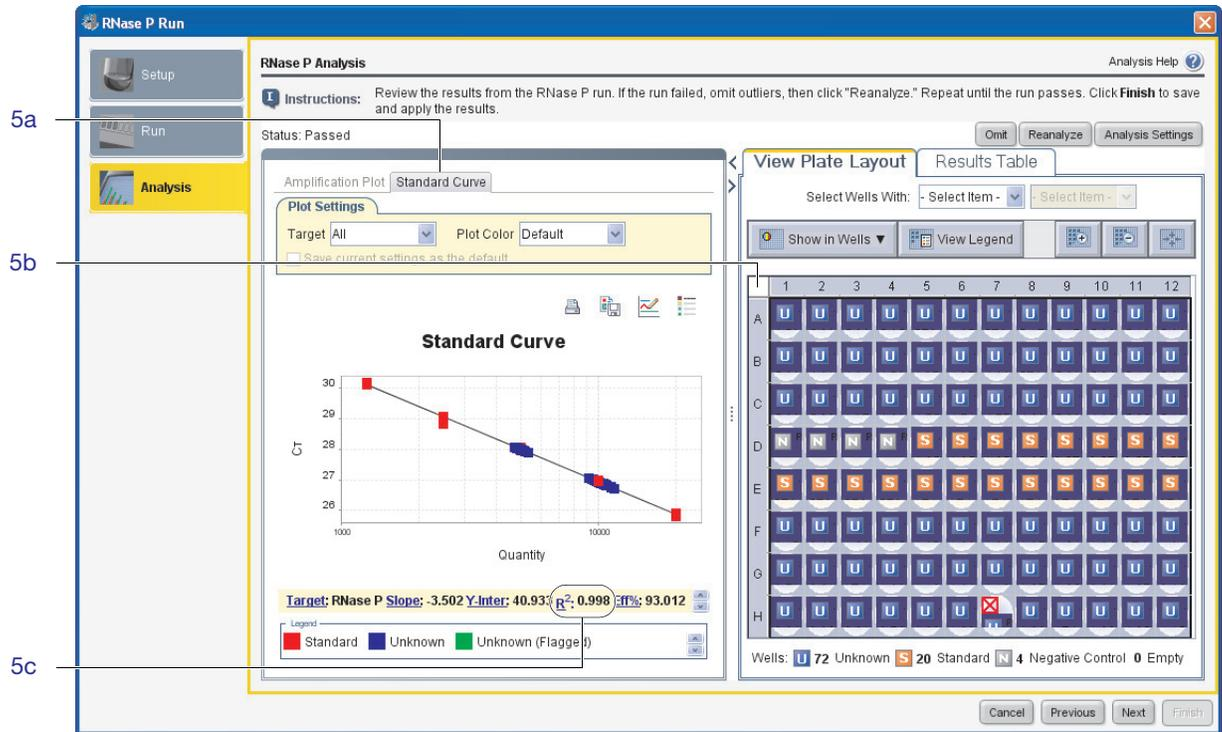
If the status of the RNase P Run is “**Failed**” after performing [steps 2](#) through [4](#), repeat the RNase P experiment using a different RNase P plate. If the problem persists, contact Applied Biosystems as explained in “[How to Obtain Support](#)” on page [x](#).

The screenshot displays the RNase P Run software interface. On the left, a sidebar shows 'Setup', 'Run', and 'Analysis' buttons. The main window is titled 'RNase P Analysis' and shows 'Status: Passed'. The 'Amplification Plot' (3b) shows a graph of Ct vs Well Number, with a single outlier well circled in red. The 'View Plate Layout' (3a) shows a 96-well plate with wells containing Unknown (U), Standard (S), and Negative Control (N) samples. A red circle highlights a well in row D, column 11, which is circled in red in the plot. The 'Results Table' (3c) shows a table with columns for 'Select Wells With:' and 'Item'. The 'Reanalyze' button (4) is visible in the top right corner. The bottom status bar shows 'Wells: U 72 Unknown S 20 Standard N 4 Negative Control 0 Empty'.

Notes _____

5. Review the standard curve:
 - a. Select the **Standard Curve** tab.
 - b. Click the upper-left corner of the Plate Layout to select all wells.
 - c. Verify that the R2 value is greater than or equal to 0.990.

If the R2 value is less than 0.990, repeat the RNase P experiment using a different RNase P plate. If the problem persists, contact Applied Biosystems as explained in “How to Obtain Support” on page x.



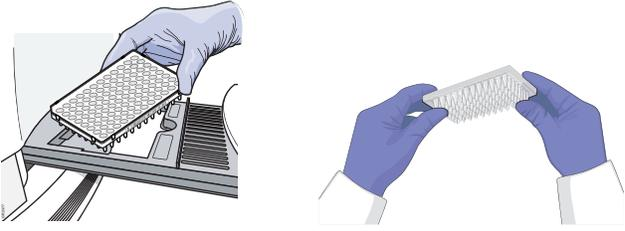
6. Click **Next**, then remove the calibration plate.

WARNING PHYSICAL INJURY HAZARD. During instrument operation, the sample block can be heated to 100 °C. Before performing the following procedure, be sure to wait until the sample block reaches room temperature.

- a. Push the tray door to open it.
 - b. Remove the calibration plate.
 - c. Push the tray door to move it into the instrument.
7. Discard the plate.
 8. Click **Finish**, then click **Yes** when prompted to save the experiment.

Notes

Troubleshoot the RNase P Experiment

Problem/Symptom	Possible Cause	Action
More than the maximum number of outliers are present in RNase P data	<ul style="list-style-type: none"> Possible contamination Pipetting inaccuracy 	Contact your Applied Biosystems service and sales representative to order a replacement TaqMan® RNase P Instrument Verification Plate. If the replacement RNase P plate fails, contact Applied Biosystems technical support or your service representative for further assistance.
The RNase P plate verification run failed	<ul style="list-style-type: none"> Insufficient centrifugation Defective plate seal 	<p> WARNING PHYSICAL INJURY HAZARD. During instrument operation, the sample block can be heated to 100 °C. Before performing the following procedure, be sure to wait until the sample block reaches room temperature.</p> <ol style="list-style-type: none"> Unload the RNase P plate from the instrument: <ol style="list-style-type: none"> Push the tray door to open it. Remove the RNase P plate from the tray. Push the tray back into the instrument. <p></p> <ol style="list-style-type: none"> Hold the plate up to a light source, and verify that all wells contain the same volume of fluid. <p>If there are differences in fluid volumes, check the heat seal of the wells with lower volumes for signs of damage or evaporation.</p> <p>Also, compare the position of the wells that have lower volumes with the outliers that you have removed from the plate. If the well positions coincide, the heat seal on the plate may be defective and resulted in the evaporation of the associated samples.</p> Contact your Applied Biosystems service and sales representative to order a replacement TaqMan® RNase P Instrument Verification Plate. If the replacement RNase P plate fails, contact Applied Biosystems technical support or your service representative for further assistance.

Notes _____

6

User-Performed Maintenance

This chapter covers:

■ Monitor the 7500/7500 Fast System	56
■ View the Instrument Log	56
■ Monitor the Instrument Status	57
■ Monitor the Lamp Status	58
■ Decontaminate the Sample Block	59
■ Monitor the 7500/7500 Fast System	58
■ Replace the Halogen Lamp	63
■ Replace the Instrument Fuses	66
■ Update the Windows Operating System	67
■ Update the 7500 Software	68

Note: For more information about any of the topics discussed in this guide, access the Help from within Applied Biosystems 7500/7500 Fast Real-Time PCR System Software v2.0 by pressing **F1**, clicking  in the toolbar, or selecting **Help ▶ 7500 Software Help**.

Notes _____

Monitor the 7500/7500 Fast System

You can monitor the state of the Applied Biosystems 7500/7500 Fast Real-Time PCR System using the Function Test, Lamp Status/Replacement, and Instrument Log tools of the 7500 software. The tools enable you to assess the health of the 7500/7500 Fast system, check the replacement status of the instrument lamp, and view a recent history of instrument activity.

View the Instrument Log

Use the Instrument Log to view the recent event history of the 7500/7500 Fast system. The log displays the major instrument activity for either the most recent 25 runs (including calibrations), or the events that pertain only to a specific EDS file.

Display the Instrument Log

- In the 7500 software, select **Instrument ▶ Instrument Events Log**.
- In the Instrument Events Log dialog box, select either:
 - System Log** – To view events that occurred during the 25 most recent runs (experiments) or calibrations.
 - Document Log** – To view events that pertain only to the experiment currently open in the 7500 software.
- If necessary, modify the data displayed by the table by filtering the data and adding or removing columns.

To...	Action
Filter the data in the events table	<ol style="list-style-type: none"> In the Filter drop-down list, select a property. Enter the appropriate conditions into the drop-down lists and fields that appear automatically. Click Apply to filter the data. <p>Note: To reset the log, select Filter ▶ Show All Records.</p>
Add or remove columns to/from the events table	Click Show Columns , then select the column desired column in the drop-down list.
Sort the data in the events table	Click the heading of the column of interest once to sort the data in ascending order. Click the column heading again to sort the data in reverse order.
Export the contents of the instrument event log	<ol style="list-style-type: none"> Select the rows of interest in the event table, Press Ctrl+C to copy the data. Paste the data into a spreadsheet application or a text file. <p>Note: The software exports the data in tab-delimited format.</p>

- When you finish viewing the events, click the close box to close the dialog box.

Notes _____

Monitor the Instrument Status

Use the Function Test dialog box to perform a high-level diagnostic of the major 7500/7500 Fast system components. In general, you need not perform the function tests unless you experience a suspected hardware failure, or you are instructed to do so by an Applied Biosystems representative.

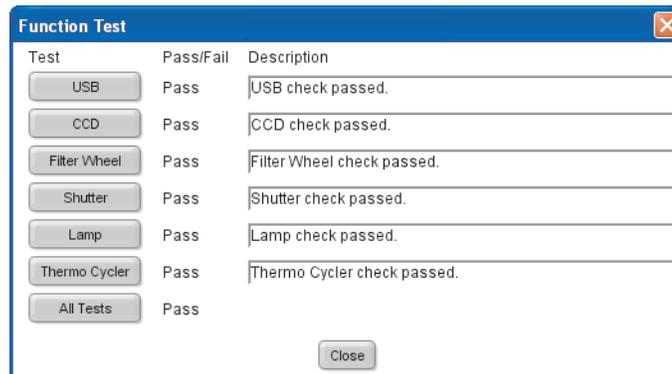
Perform Function Tests of System Components

1. In the 7500 software, select **Instrument ▶ Function-Test**.
2. Perform function tests as needed.

To test:

- All system components – Click **All Tests**, then wait for the software to perform all of the function tests.
- One or more specific components – Click one or more of the following.
 - USB** – Tests the universal serial bus (USB) connection between the 7500/7500 Fast instrument and computer. The test passes if the 7500 software can establish communication with the 7500/7500 Fast instrument.
 - CCD** – Tests the CCD camera in the 7500/7500 Fast instrument. The test passes if the camera can capture an image.
 - Filter Wheel** – Tests the filter wheel in the 7500/7500 Fast instrument. The test passes if the filter wheel controller is running the correct version of firmware.
 - Shutter** – Tests the optic shutter of the 7500/7500 Fast instrument. The test passes if the shutter controller is running the correct version of firmware.
 - Lamp** – Tests the halogen lamp of the 7500/7500 Fast instrument. The test passes if the lamp controller is running the correct version of firmware.
 - Thermal Cycler** – Tests the thermal cycler sample block in the 7500/7500 Fast instrument. The test passes if the thermal cycler controller is running the correct version of firmware.

When the 7500 software completes a test, the software reports the pass/fail status of the test and provides a description of the outcome.



3. When you finish testing, click **Close**.

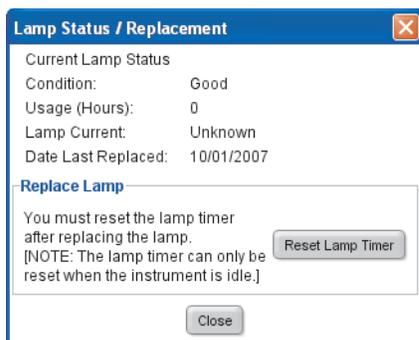
Notes

Monitor the Lamp Status

Use the Lamp Status/Replacement dialog box to monitor the status of the halogen lamp that the 7500/7500 Fast instrument uses to illuminate samples during runs.

Check the Lamp Status

In the 7500 software, select **Instrument** ▶ **Lamp Status/Replacement** to determine the status of the halogen lamp.



The Lamp Status/Replacement dialog box displays:

- **Condition** – Indicates one of the following conditions:
 - **Good** – The lamp is functioning well and does not need to be replaced. Click **Close**.
 - **Failed** – The lamp bulb must be replaced. Click **Close**, then replace the lamp as explained in “[Replace the Halogen Lamp](#)” on page 63.
 - **Change Soon** – The lamp usage is above 2000 hrs; Applied Biosystems recommends that you replace it. Click **Close**, then decide whether or not to replace the lamp. If you choose to replace the lamp, see “[Replace the Halogen Lamp](#)” on page 63.
- **Usage (Hours)** – The total number of hours that the lamp has been illuminated.
- **Lamp Current** – The output current of the lamp in amperes (A). Low current can indicate a potential future failure of the lamp.
- **Date Last Replaced** – The date of the last lamp replacement.

Warnings The 7500 software can display the following warnings before or during a run:

Message	Description
Warning – Cannot detect sufficient current from lamp. Either lamp is not installed properly or needs to be replaced.	The lamp current is below the acceptable level at the start of the run. You cannot proceed with the run until you replace the halogen bulb as explained in “ Replace the Halogen Lamp ” on page 63.
Warning – Cannot detect sufficient current from lamp. Either lamp is not installed properly or needs to be replaced.	The 7500 software stopped the run because the lamp current decreased below the acceptable level during the run. You cannot proceed with the run until you replace the halogen bulb as explained in “ Replace the Halogen Lamp ” on page 63. Click OK in the message box, inspect the Instrument Log, then replace the lamp bulb.
Warning - The lamp usage has exceeded 2000 hr. We recommend replacing the lamp soon to ensure optimal assay performance.	The lamp usage exceeds 2000 hr at the start of a run. Click Cancel Run , then replace the lamp, or click Continue Run . Rerun ROI, Background, Optical and Dye calibrations.

Notes

Decontaminate the Sample Block

Perform the following procedure to eliminate fluorescent contaminants from the sample block of the 7500/7500 Fast instrument. Fluorescent contamination is generally evident in failed background runs where one or more wells consistently exhibit abnormally high signals.

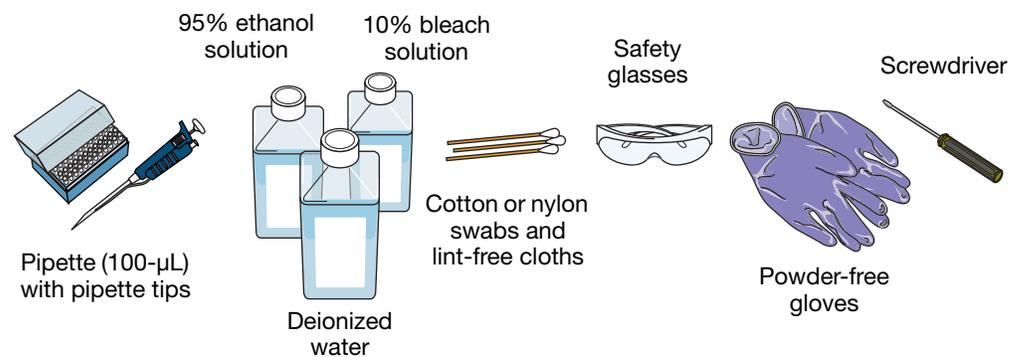
CAUTION **PHYSICAL INJURY HAZARD.** Do not remove the instrument cover. There are no components inside the 7500/7500 Fast system that you can safely service yourself. If you suspect a problem, contact an Applied Biosystems Service Representative.

CAUTION **PHYSICAL INJURY HAZARD.** During instrument operation, the sample block can be heated to 100 °C. Before performing the following procedure, be sure to wait until the sample block reaches room temperature.

CAUTION Before using a cleaning or decontamination method other than those recommended by the Applied Biosystems, verify with Applied Biosystems that the proposed method will not damage the equipment.

Time Required 30 min

Materials Required



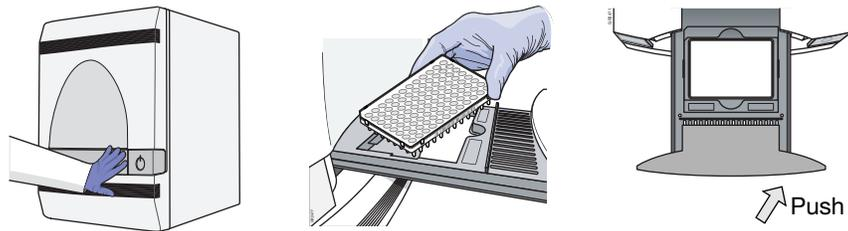
Clean the Sample Block

IMPORTANT! Wear powder-free gloves when you perform this procedure.

1. Identify the contaminated wells of the sample block (see “[How to Identify Contamination](#)” on page 30).
2. Remove the plate and the tray holder from the 7500/7500 Fast instrument:
 - a. Push the tray door to open it.
 - b. Remove the plate and the tray holder.

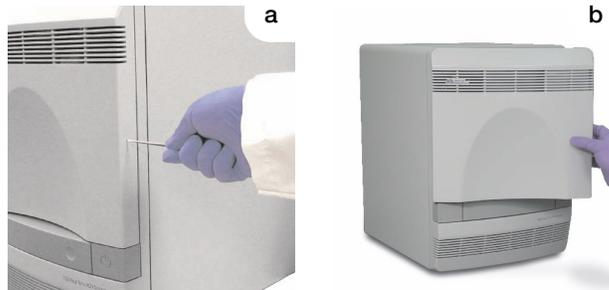
Notes _____

- c. Close the tray door. Apply pressure to the right side of the tray door at an angle.



3. Manually raise the block:

- a. In the 7500 software, select **Instrument ▶ Instrument Maintenance Manager**.
 - b. In the ROI tab of the Instrument Maintenance Manager, click **Start Manual Calibration**.
 - c. In the ROI Inspector dialog box, click **Move Block**.
 - d. When the ROI Inspector dialog box displays “Block Down,” click **Done**.
4. Power off, then unplug the 7500/7500 Fast instrument. Allow it to cool for 15 min.
5. Open the access door to the 7500/7500 Fast instrument.
- a. Insert a thin screwdriver into the keyhole on the edge of the access door, then push to unlatch the door.
 - b. Open the access door.



6. Lift the latch, then push the heated cover door to the back of the instrument.

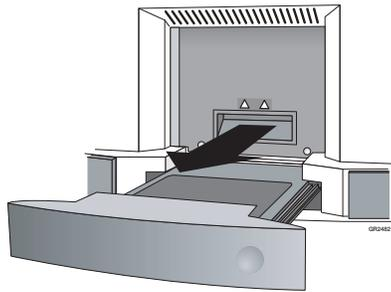


Notes _____

7. Clean the contaminated wells of the sample block using a small volume of deionized water:
 - a. Pipette a small volume of deionized water into each contaminated well.
 - b. Pipette the water up and down several times to rinse the well.
 - c. Pipette the water to a waste beaker.
 - d. Using a cotton swab, scrub inside of each contaminated well.
 - e. Using a lint-free cloth, absorb the excess deionized water.



8. Pull the heated cover door to the front of the 7500/7500 Fast instrument. Lift the latch, then secure the heated cover door to the cross bar.



9. Close the access door to the 7500/7500 Fast instrument.



10. Plug in, then power on the 7500/7500 Fast system.
11. Verify that you have eliminated the contamination by performing a background calibration run (see [“Perform the Background Calibration”](#) on page 20).

Notes _____

12. If the contamination remains, repeat [steps 1 through 6](#), then clean the contaminated wells of the sample block using 95% ethanol solution:
 - a. Pipette a small volume of 95% ethanol solution into each contaminated well.
 - b. In each contaminated well, pipette the solution up and down several times to rinse the well.
 - c. Pipette the ethanol solution to a waste beaker.

IMPORTANT! Always use deionized water to rinse wells after cleaning with bleach or ethanol solution.

13. Repeat [steps 7 through 11](#) to rinse the wells of the sample block and to verify that you have eliminated the contamination.
14. If the contamination remains, repeat [steps 1 through 6](#), then clean the contaminated wells of the sample block using 10% bleach solution:
 - a. Pipette a small volume of 10% bleach solution into each contaminated well.
 - b. In each contaminated well, pipette the solution up and down several times to rinse the well.
 - c. Pipette the bleach solution to a waste beaker.

IMPORTANT! Always use deionized water to rinse wells after cleaning with bleach or ethanol solution.

15. Repeat [steps 7 through 11](#) to rinse the wells of the sample block and to verify that you have eliminated the contamination.
If the contamination remains, contact Applied Biosystems support (see [“How to Obtain Support”](#) on page x).
16. Ensure that the heated cover door is completely closed and latched. If it is not, the 7500 software displays an error message.

Notes _____

Replace the Halogen Lamp

Replace the halogen lamp after approximately 2000 hr of life.

WARNING **PHYSICAL INJURY HAZARD.** The 7500/7500 Fast system and lamp are hot! The lamp can become very hot while in use. Allow the lamp to cool for 15 min and put on protective, powder-free gloves before handling it.

CAUTION **PHYSICAL INJURY HAZARD.** Wear disposable, powder-free gloves when handling the lamp to prevent burns and to prevent shortening the life of the replacement lamp.

WARNING **WARNING.** This instrument is designed for 12 V, 75 W halogen lamps only.

Time Required 30 min

Materials Required



Halogen bulb
(12 V, 75 W)



Screwdriver,
small



Powder-free
gloves



Safety
glasses

Replace the Lamp

IMPORTANT! Wear powder-free gloves when you handle the lamp.

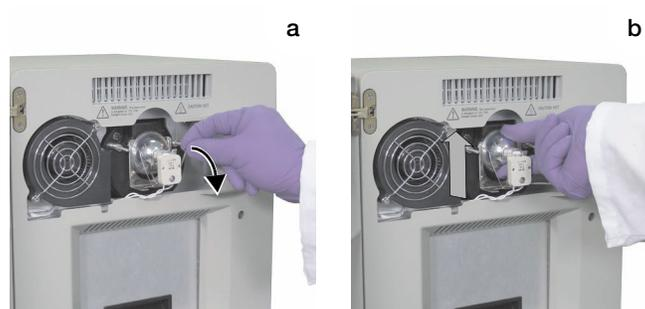
1. Power off, then unplug the 7500/7500 Fast system. Allow the instrument to cool for 15 min.
2. Open the access door to the 7500/7500 Fast system:
 - a. Insert a thin screwdriver into the keyhole on the edge of the access door, then push to unlatch the door.
 - b. Open the access door.



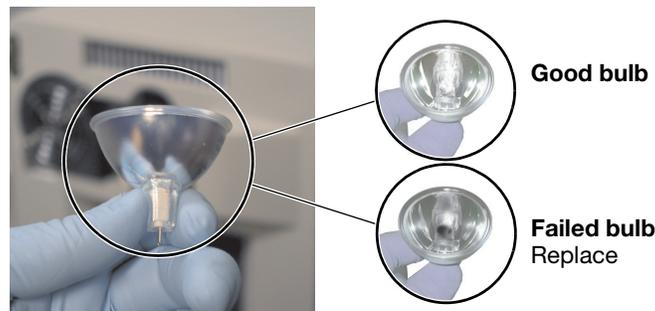
Notes

3. Remove the lamp from the instrument:
 - a. Slide the lamp release lever downward.
 - b. Firmly grasp the lamp and lift it up and out of the slotted mount.

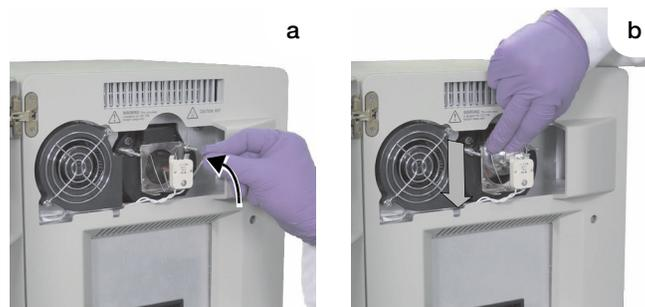
IMPORTANT! Do not touch the lamp without powder-free gloves. Finger prints shorten the lamp life.



4. Inspect the lamp for signs of failure (carbon typically coats the inside of failed lamps).



5. Install the new lamp into the instrument:
 - a. Slide the lamp release lever upward.
 - b. Firmly grasp the lamp, place it into the slotted mount, then carefully slide the lamp downward into place.

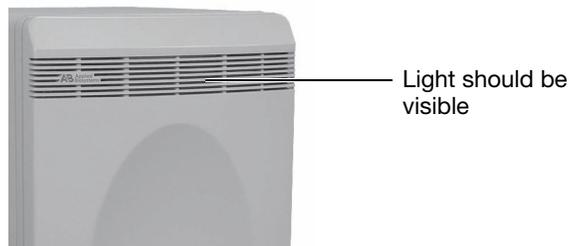


Notes _____

6. Close the access door.



7. Plug in and power on the 7500/7500 Fast system.
8. Open the ROI Inspector dialog box:
 - a. In the 7500 software, select **Instrument** ▶ **Instrument Maintenance Manager**.
 - b. In the ROI tab of the Instrument Maintenance Manager, click **Start Manual Calibration**.
9. In the ROI Inspector dialog box, select **Lamp Control** ▶ **Idle**.
10. While the instrument is running, look through grating of the access door and verify that the lamp is illuminated, then click **Done**.



11. If the lamp is illuminated, select **Instrument** ▶ **Lamp Status/Replacement** in the 7500 software, click **Reset Lamp Timer**, then click **OK**.

If the lamp is not illuminated, the replacement halogen lamp may be defective. Replace the lamp again. If the second lamp does not illuminate, check the instrument fuses for failure (see [page 66](#)).
12. Perform the following calibrations after replacing the lamp. See:
 - [Chapter 2, Perform the Regions of Interest \(ROI\) Calibration](#)
 - [Chapter 3, Perform the Background Calibration and Optical Calibration](#)
 - [Chapter 4, Perform the Dye Calibration](#)
 - [Chapter 5, Verify the Instrument Performance](#)

Notes _____

Replace the Instrument Fuses

Replace the 7500/7500 Fast instrument fuses when the fuses fail.

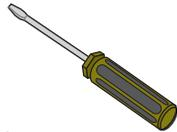
 **CAUTION FIRE HAZARD.** For continued protection against the risk of fire, replace fuses only with listed and certified fuses of the same type and rating as those currently in the instrument.

Time Required 30 min

Materials Required



Fuses (2),
12.5 A, 250 V,
5 x 20 mm



Flathead
screwdriver



Powder-free
gloves



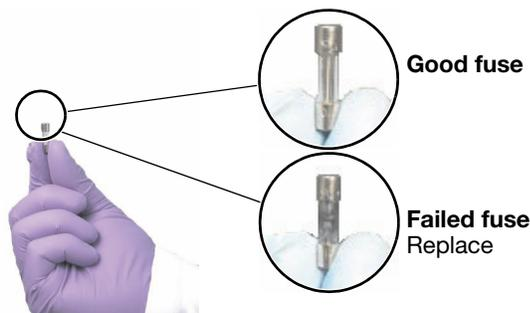
Safety
glasses

Replace the Fuses

1. Power off the instrument, then unplug it.
2. Using a flat-head screwdriver, unscrew and remove the fuse holders from the instrument.



3. Remove each fuse from its fuse holder and inspect it for damage. Carbon typically coats the inside of failed fuses.



Notes _____

4. Replace each failed fuse with a 12.5 A, 250 V, 5 × 20-mm fuse.

Note: The voltage and amperage ratings are on the fuse holder.

5. Install the fuse holder.



6. Plug in, then power on the instrument.

The installation is successful if the instrument powers on.

Note: Fuse failure can result from fluctuations in the supplied power to the instrument. To prevent further failures, consider installing an electrical protective device, such as a UPS or other surge protector. For more information about fuses, see the *7500/7500 Fast Site Preparation Guide*.

Update the Windows Operating System

Do not upgrade or update the Microsoft Windows® operating system of the computer running the 7500 software without first consulting the software release notes or the Applied Biosystems website. Future versions of the Windows® operating system and updates to the operating system can conflict with the 7500 software.

Determine Compatibility of an Upgrade or Update

1. Open **D:\Applied Biosystems\7500 Software v2.0**, double-click **release-notes.html**, then read the *7500 Software Release Notes* for the compatibility of interest.
2. If the release notes do not mention the compatibility, use an internet browser to visit **www.appliedbiosystems.com**, then search the website for the compatibility of interest.
3. If the website does not contain the information of interest, contact Applied Biosystems (see “[How to Obtain Support](#)” on page x).

Notes _____

Update the 7500 Software

If you want to update the 7500 software, prepare your computer by exporting the application libraries and backing up your experiment files.

Visit the Applied Biosystems Website

You can obtain 7500 software updates directly from the service section of the Applied Biosystems website. For the latest services and support information for the 7500/7500 Fast instrument:

1. Go to <https://www2.appliedbiosystems.com/support/software/>
2. In the Software Downloads page, select the appropriate instrument from the drop-down list.
3. In the Software Downloads page for your instrument, click **Updates - Patches**.

The website opens the page describing the latest software updates for the 7500 software.

Prepare for the Upgrade

Before updating the 7500 software:

1. Back up the application libraries:
 - a. In the main menu of the 7500 software, select **Tools** ▶ *<desired library>*.
 - b. When the library dialog box opens, select the element(s) that you want to export, then click **Export**.
 - c. In the Export dialog box, click **Save** to archive the selected records.
 - d. Repeat [steps 1a](#) through [1c](#) for the remaining libraries to archive them.
2. Back up all experiment files by creating a copy of the directory that you are using to store files.

The default directory for experiments is:

D:\Applied Biosystems\7500\experiments

Notes _____



Store, Move, and Install the 7500/7500 Fast System

A

This chapter covers:

- Store the 7500/7500 Fast System 70
- Move the 7500/7500 Fast System 71
- Set Up the 7500/7500 Fast System 73

Note: For more information about any of the topics discussed in this guide, access the Help from within Applied Biosystems 7500/7500 Fast Real-Time PCR System Software v2.0 by pressing **F1**, clicking  in the toolbar, or selecting **Help ▶ 7500 Software Help**.

Notes

Store the 7500/7500 Fast System

The Applied Biosystems 7500/7500 Fast Real-Time PCR System can be powered off and stored for extended periods of time. The length of the period of inactivity determines the method you use to power off the instrument.

Time Required 5 min

Materials Required ABI PRISM® Optical 96-Well Reaction Plate or Optical 96-Well Fast Plate (unused)

Prepare the Instrument

1. Open the instrument tray door.
2. If the tray contains a plate, remove it.
3. If you plan to store the 7500/7500 Fast instrument for more than a week or you plan to move the instrument, load an unused plate into the tray.

Note: The empty plate protects the internal components of the 7500/7500 Fast instrument during transport or during periods of inactivity lasting more than a week.

4. Push the tray door to move it into the instrument.



5. Press the instrument power button.
6. Power off the computer and monitor:
 - a. Select **Start ▶ Shut Down**.
 - b. In the Shut Down Windows dialog box, select **Shut Down**.
 - c. Power off the monitor.

Notes _____

Move the 7500/7500 Fast System

Perform this procedure to safely move the 7500/7500 Fast instrument short distances (for example, between laboratories of the same building).



CAUTION

PHYSICAL INJURY HAZARD. Do not attempt to lift the instrument or any other heavy objects unless you have received related training. Incorrect lifting can cause painful and sometimes permanent back injury. Use proper lifting techniques when lifting or moving the instrument. At least 2 people are required to lift the 7500/7500 Fast system.

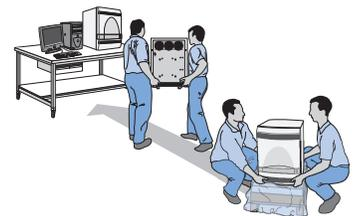
IMPORTANT! Moving your Applied Biosystems 7500/7500 Fast Real-Time PCR System can create subtle changes in the alignment of the instrument optics.

Materials Required

ABI PRISM® Optical 96-Well Reaction Plate and Optical 96-Well Fast Plate (unused)

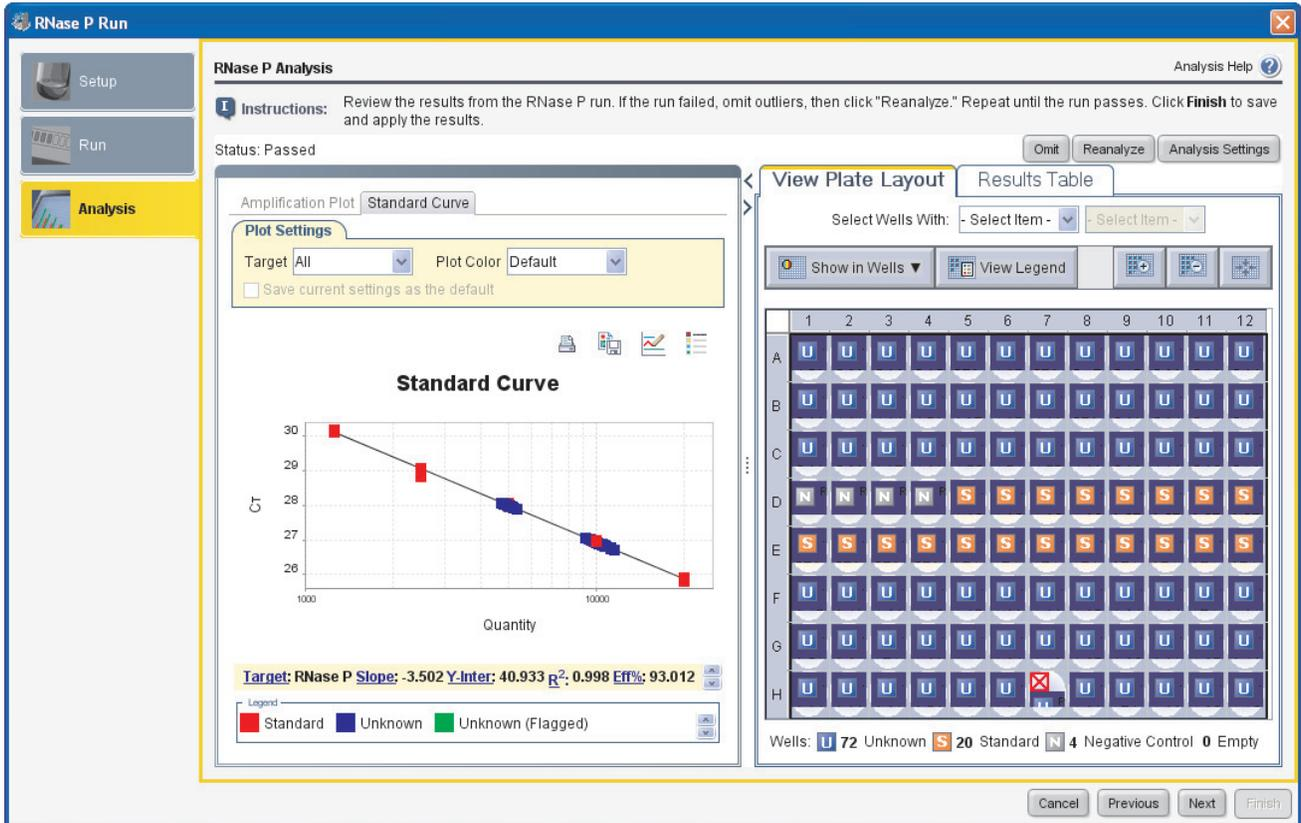
Prepare the Instrument

1. Load the empty reaction plate into the 7500/7500 Fast instrument.
2. Using the ROI Inspector, manually raise the sample block:
 - a. In the 7500 software, select **Instrument ▶ Instrument Maintenance Manager**.
 - b. In the ROI tab of the Instrument Maintenance Manager, click **Start Manual Calibration**.
 - c. In the ROI Inspector dialog box, click **Block Up**.
3. Power off the 7500/7500 Fast instrument and computer.
4. Disconnect all 7500/7500 Fast system components.
5. Move the 7500/7500 Fast system according to the following guidelines:
 - Verify that the surface on which you will place the instrument can support at least 54.5 kg (120 lbs).
 - Verify that the pathway to the final position of the instrument is clear of obstructions.
 - Keep your spine in a good neutral position.
 - Bend at the knees and lift with your legs.
 - Do not lift an object and twist your torso at the same time.
 - Coordinate your intentions with your assistant before lifting and carrying.



Notes

6. Reconnect the components of the 7500/7500 Fast system (see “Set Up the 7500/7500 Fast System” on page 73).
7. Run a TaqMan® RNase P Instrument Verification Plate (see page 45).
 - a. If the run passes, recalibrations are not necessary.
 - b. If the run fails, perform steps 8 through 12 to recalibrate the instrument.



8. Perform an ROI calibration (see page 13).
9. Perform a background calibration (see page 23).
10. Perform an optical calibration (see page 25).
11. Perform a dye calibration (see page 32).
12. Perform an instrument verification run (see page 45).

Notes _____

Set Up the 7500/7500 Fast System

Set Up the Computer

Refer to the *Applied Biosystems Real-Time System Computer Setup Guide* for information on setting up a computer for use with the 7500/7500 Fast instrument.

Set Up the 7500/7500 Fast Instrument

IMPORTANT! Do not connect the USB cable to the 7500/7500 Fast instrument until you are instructed to do so by this guide.

Materials Required

- Phillips screwdriver (small and thin)
- Power cord

Set Up the 7500/7500 Fast System

1. Prepare the installation site as described in the *Applied Biosystems 7500/7500 Fast Real-Time PCR System Site Preparation Guide*.
2. Open the access door to the 7500/7500 Fast instrument.
 - a. Insert a thin screwdriver into the keyhole on the edge of the access door, then push to unlatch the door.
 - b. Open the access door.
3. Verify that the heated cover assembly is pulled fully toward the front of the instrument. If the 7500/7500 Fast system has a heated cover latch installed, check that the latch is in a locked position.
4. Inspect the instrument for damage caused by the transportation of the 7500/7500 Fast system.

If the instrument is damaged, record the location and appearance of the damage, then contact Applied Biosystems technical support or your service representative for assistance.
5. Close the access door.
6. Connect the power cord to the 7500/7500 Fast instrument, then to the wall receptacle.

Note: Power cords for different voltages are provided in the packing kit. Connect the cord with the receptacle appropriate for your voltage, then discard the remaining power cords.

7. Press the power button at the lower right front panel, then wait for the 7500/7500 Fast instrument to start up (about 30 sec).

Notes



8. When the Power status light on the lower left front panel is lit, push the tray door to open it.
9. Remove the packaging plate from the tray and set it aside.
10. Close the tray door, then press the power button again to power off the instrument.

Note: Install any additional hardware.

IMPORTANT! Do not connect the USB cable to the 7500/7500 Fast instrument at this time.

11. Verify that the 7500 software is installed to the computer.
If the computer does not have the 7500 software, use the Applied Biosystems 7500/7500 Fast Real-Time PCR System Software v2.0 Software CD to install the software.
12. Once you verify that the computer contains the 7500 software, connect the USB cable to the 7500/7500 Fast instrument.

Notes _____

B

Create a Custom Dye Plate

The Applied Biosystems 7500/7500 Fast Real-Time PCR System can be used to run assays designed with custom dyes (dyes not manufactured by Applied Biosystems). Custom dyes must fluoresce within the 520 to 650 nm spectral range measured by the 7500/7500 Fast instrument.

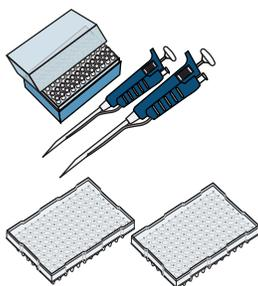
Before Using Custom Dyes

Before using custom dyes with the 7500/7500 Fast instrument, you must:

- Determine optimum dye concentration
- Create a custom dye plate
- Add the custom dye to the software
- Perform a dye calibration (see [Chapter 4](#), “Perform the Dye Calibration.”)

Materials Required

Pipettors and pipette tips (200- μ L and 1000- μ L)

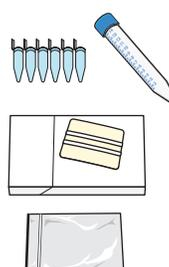


ABI PRISM® Optical 96-Well Reaction Plate and Optical 96-Well Fast Plate

Water, deionized



Tubes (2-mL and 10-mL)



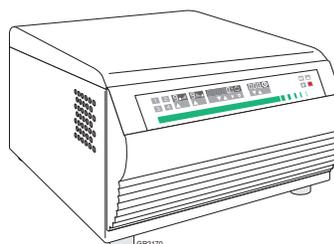
Optical Adhesive Cover or Optical Flat Caps

Safety glasses



Powder-free gloves

Custom dye(s)



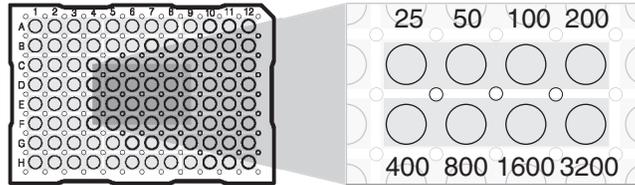
Centrifuge with plate adapter

Notes

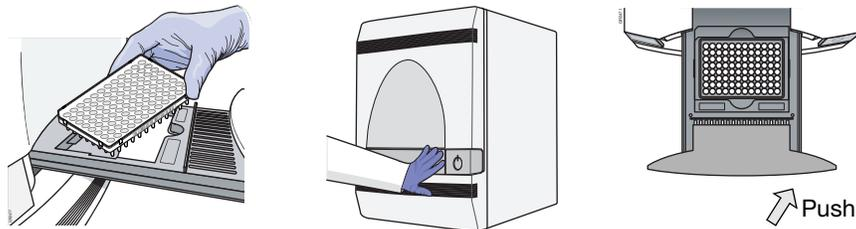
Determine Optimum Dye Concentration

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

1. In the center wells of a 96-well plate, prepare a dilution series of the custom dye (for example, 25, 50, 100, 200, 400, 800, 1600, and 3200 nM) using 50- μ L volumes for the 7500 system or 20- μ L volumes for the 7500 Fast system.



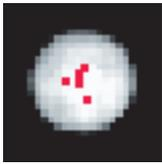
2. Seal the wells of the reaction plate using an optical adhesive cover.
3. Load the prepared reaction plate:
 - a. Push the tray door to open it.
 - b. Load the dye plate into the plate holder.
 - c. Push the tray door to close it.



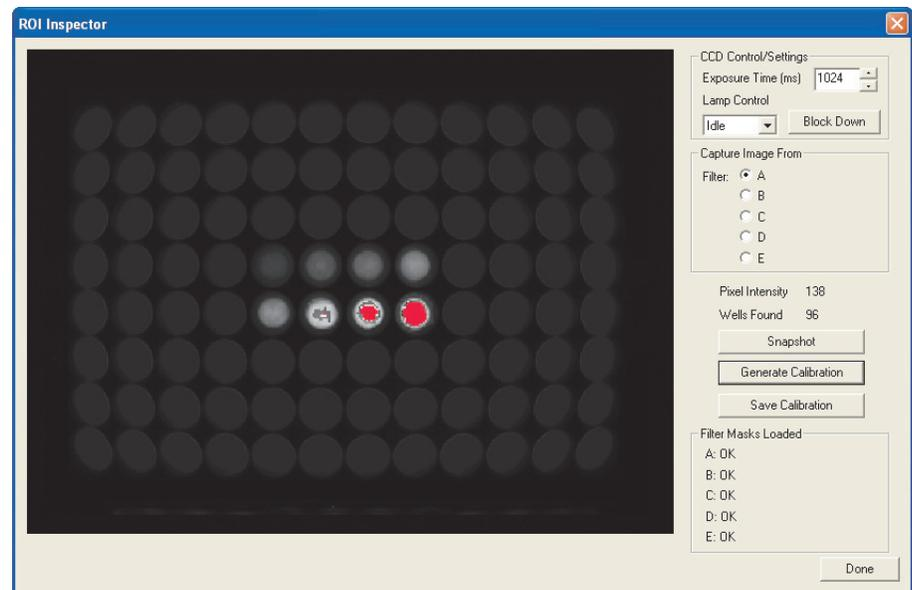
4. In the 7500 software, select **Instrument** ► **Instrument Maintenance Manager**.
5. In the ROI tab of the Instrument Maintenance Manager, click **Start Manual Calibration**.
6. In the ROI Inspector, create the ROI image for each filter, beginning with Filter 1:
 - a. In the Exposure Time field, enter **1024**.
 - b. Click **Block Up**.
 - c. Select **Filter 1**.
 - d. Click **Snapshot**.

Notes _____

- e. Inspect the image for saturation.

No Saturation	Normal Saturation	Over Saturation
		
Acceptable	Not Acceptable	

- f. Record the coordinate of the well that displays the brightest possible signal without saturation. This well contains the optimal concentration of the custom dye for Filter 1.



7. Repeat steps 6c through 6f for the remaining filters.
8. After you determine the optimum concentration for each filter, determine the optimum concentration for the custom dye:
 - a. Compare the results from all filters.
 - b. Select the concentration that yields the highest possible signal in all filters, but does not saturate.

Notes _____

Unload the Plate



CAUTION **PHYSICAL INJURY HAZARD.** During instrument operation, the sample block can be heated to 100 °C. Before performing the following procedure, be sure to wait until the sample block reaches room temperature.

1. In the ROI Inspector, click **Block Down**.
2. Remove the plate from the instrument:
 - a. Push the tray door to open it.
 - b. Remove the plate.
 - c. Push the tray door to move it into the instrument.

Note: If you cannot open the tray, the sample block may be in its raised position, locking the tray position. To lower the block, click **Move Down**, then click **Block Down**.

3. Click **Done**.

Create a Custom Dye Plate

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

1. Prepare 5 mL (7500 system) or 2 mL (7500 Fast system) of the custom dye at the concentration determined in [step 8 on page 77](#).
2. Pipette 50 µL (7500 system) or 20 µL (7500 Fast system) of the diluted custom dye to all wells of an optical reaction plate.
3. Seal the wells of the reaction plate using an optical adhesive cover.

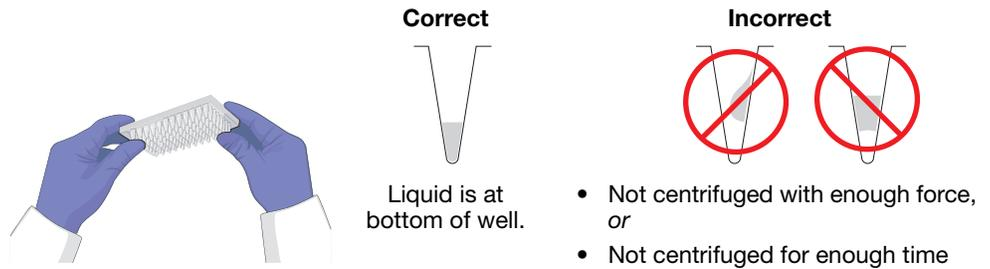


4. Centrifuge the plate for 2 min at less than 1500 rpm.

IMPORTANT! The custom dye calibration plate must be well mixed and centrifuged.

Notes _____

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



Add the Custom Dye to the Software

1. In the main screen of the 7500 software, select **Instrument ▶ Instrument Maintenance Manager**.
2. In the Instrument Maintenance Manager:
 - a. In the navigation pane, click **Dye**.
 - b. In the Dye screen, select **Custom Dye Calibration**.
 - c. Click **Start Calibration**.
3. In the Setup screen of the Dye Calibration dialog box, select a custom dye from the list or create the custom dye as follows:
 - a. Click **New Dye**.
 - b. In the Dye Manager dialog box, click **New**.
 - c. Complete the New Dye dialog box, then click **OK**.

Field/Option	Action
Name	Enter a name for the custom dye.
Wavelength	Enter the wavelength at which the dye fluoresces.
Type	Select: <ul style="list-style-type: none"> • Reporter if the dye works in conjunction with a quencher dye to report an increase of PCR product. • Quencher if the dye suppresses the fluorescence of a reporter dye until amplification of PCR product. • Both if the dye reports an increase of PCR product without the aid of a quencher dye.

- d. Click **Close**.

Notes _____

4. In the Setup screen of the Dye Calibration dialog box, enter a temperature setting for the calibration.

Note: Set the temperature to match the temperature at which you intend to collect data. For example, the temperature for all Applied Biosystems system dyes is 60 °C because data collection for TaqMan® reagents occurs during the 60 °C extension step of the PCR.

5. Select **The custom dye plate is loaded in the instrument**, then click **Next**.
6. In the Run screen, click **Start Run**, then wait for the 7500/7500 Fast instrument to complete the dye calibration.

Note: If the 7500 software displays messages during the run, troubleshoot the errors as described in [“Troubleshoot the Dye Calibration” on page 43](#).

7. When the 7500/7500 Fast instrument displays the Main Menu, unload the custom calibration plate.
8. Analyze the custom spectral calibration as explained in [“Analyze the Calibration Data” on page 39](#).

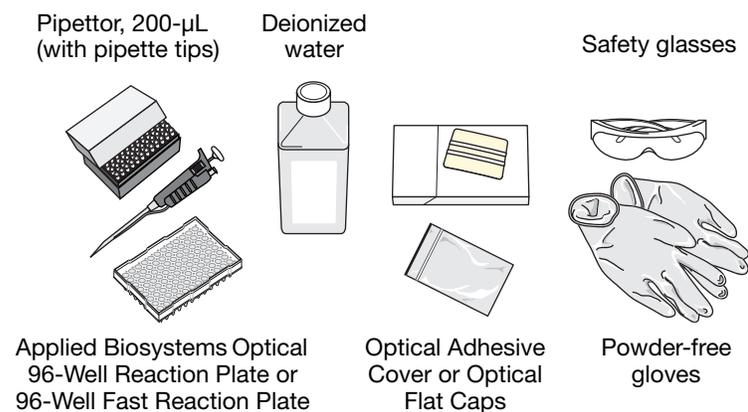
Notes _____



Create a Background Plate

Whenever possible, use a background plate that is included with the spectral calibration kit. The plates supplied in the kit contain a buffer that accurately simulates the reagents used for PCR, and, therefore, produces high-quality calibration data. However, if a background plate from a spectral calibration kit is not available, you can create one as described below.

Materials Required



Create a Background Plate

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

1. Remove an Applied Biosystems 96-Well Optical Reaction Plate or 96-Well Fast Reaction Plate from its box and place it on a clean, dry surface.
2. Aliquot 50 μ L (7500 system) or 20 μ L (7500 Fast system) of deionized water to each well of the reaction plate.
3. Seal the plate using an optical adhesive cover or optical flat caps.

Use the plate for background calibration in the same way you use a background plate from the spectral calibration kit. See [Chapter 3, “Perform the Background Calibration and Optical Calibration,”](#)

Notes



Notes _____

A

- access door, open 63
- analyze
 - background calibration 23
 - dye calibration 39–42
 - optical calibration 28
 - RNase P experiment 51–53
 - ROI calibration (automated) 12
 - ROI calibration (manual) 13–15
- Applied Biosystems
 - contacting x
 - Technical Support x
- assumptions for using this guide vii

B

- background calibration 20–24
 - analyze the data 23
 - materials required 20
 - perform 23
 - preparation 21
 - troubleshoot 30
- background fluorescence 20
- background plate
 - how to create 81
 - prepare 21
- biohazardous waste, handling xix
- biological hazard guidelines xxi

C

- calibration
 - background 20–24
 - dye 32–42
 - dye (custom) 75–80
 - optical 25–29
 - ROI 8–17
- CAUTION, description xii
- CCD function test 57
- chemical safety xvii, xviii
- chemical waste safety xix
- conventions
 - for describing menu commands vii
 - IMPORTANT! vii
 - Notes vii
 - safety xii
 - text vii
 - user attention words vii
- create

- background plate 81
 - custom dye plate 75–80
- custom dye
 - add to the software 79
 - calibration 79–80
 - plate, how to create 75–80
- CY3® dye 33, 41
- CY5® dye 33, 42

D

- DANGER, description xii
- decontamination
 - clean the sample block 59–62
 - identify contaminants 30
 - materials required 59
- documentation, related viii
- dye calibration
 - analyze the data 39–42
 - materials required 32
 - perform 36
 - preparation 35
 - spectra evaluation 34
 - spectra, 7500 41
 - troubleshoot 43

E

- electrical safety xx
- electromagnetic compatibility standards.
 - See EMC standards
- EMC standards xxii
- ergonomics, safety xxi

F

- FAM™ dye 33, 41, 46
- Fast plate
 - description 9
 - do not vortex 9
- Filter Wheel function test 57
- fluorescence, background 20
- function tests 57
- fuse replacement 66–67
 - materials required 66

G

- guidelines
 - chemical safety xviii

chemical waste disposal [xix](#)
 chemical waste safety [xix](#)

H

halogen lamp. *See* lamp
 hazard icons. *See* safety symbols, on instruments
 hazard symbols. *See* safety symbols, on instruments
 hazards. *See* safety
 Help system, accessing [ix](#)

I

IMPORTANT, description [xii](#)
 installation
 halogen lamp [63–65](#)
 instrument fuses [66–67](#)
 operating system updates [67](#)
 specification [47](#)
 installation category [xx](#)
 instrument
 background calibration [20–24](#)
 dye calibration [32–42](#)
 dye calibration (custom) [75–80](#)
 load [49](#)
 move [71](#)
 optical calibration [25–29](#)
 prepare for storage [70](#)
 RNase P experiment [46–53](#)
 ROI calibration [8–17](#)
 setup [73](#)
 instrument fuses
 replacement [66–67](#)
 specification [66](#)
 instrument log, view [56](#)
 instrument operation, safety [xvi](#)

J

JOE™ dye [33, 41](#)

L

Lamp function test [57](#)
 lamp replacement [63–65](#)
 failed lamp, example [64](#)
 lamp removal [64](#)
 materials required [63](#)
 safety [63](#)
 troubleshoot [65](#)
 lamp status, view [58](#)
 lamp timer, reset [58](#)
 lamp, life span [63](#)
 load, instrument [49](#)

M

materials required

background calibration [20](#)
 decontamination [59](#)
 dye calibration [32](#)
 fuse replacement [66](#)
 lamp replacement [63](#)
 optical calibration [25](#)
 RNase P experiment [46](#)
 ROI calibration [8](#)
 move, instrument [71](#)
 moving and lifting, safety [xvi](#)
 moving parts, safety [xx](#)
 MSDSs
 description [xvii](#)
 obtaining [x, xvii](#)

N

NED™ dye [33, 41](#)

O

omit well [51](#)
 online Help. *See* Help system
 operating system, update [67](#)
 optical calibration [25–29](#)
 analyze the data [28](#)
 materials required [25](#)
 perform [28](#)
 preparation [25](#)
 troubleshoot [30](#)
 outliers [51](#)
 overvoltage category (rating) [xx](#)

P

perform
 background calibration [23](#)
 dye calibration [36](#)
 optical calibration [28](#)
 RNase P experiment [49](#)
 ROI calibration [12](#)
 physical hazard safety [xx](#)
 plate
 background [21](#)
 dye [35](#)
 Fast [9](#)
 RNase P [46, 48](#)
 ROI [9, 25](#)
 standard [9](#)
 preparation
 background calibration [21](#)
 dye calibration [35](#)
 optical calibration [25](#)
 RNase P experiment [48](#)
 ROI calibration [9](#)
 ROI plate [25](#)

R

- R2 value 53
- radioactive waste, handling xix
- remove outliers from RNase P runs 51
- repetitive motion, safety xxi
- replace
 - halogen lamp 63–65
 - instrument fuses 66–67
- reset lamp timer 58
- RNase P experiment 46–53
 - analyze 51–53
 - materials required 46
 - perform 49
 - preparation 48
 - prepare 48
 - R2 value 53
 - troubleshoot 54
- ROI calibration 8–17
 - analyze the data (automated) 12
 - analyze the data (manual) 13–15
 - materials required 8
 - perform 12
 - preparation 9
 - saturation examples 13
 - successful 14
 - troubleshoot 17
- ROX™ dye 33, 42

S

- safety
 - before operating the instrument xvi
 - biological hazards xxi
 - chemical xvii
 - chemical waste xix
 - conventions xii
 - electrical xx
 - ergonomic xxi
 - guidelines xviii, xix
 - instrument operation xvi
 - moving and lifting instrument xvi
 - moving parts xx
 - moving/lifting xvi
 - physical hazard xx
 - repetitive motion xxi
 - standards xxii
 - workstation xxi
- safety labels, on instruments xiv
- safety standards xxii
- safety symbols, on instruments xiii
- saturation, ROI calibration 13
- service pack, updates 67
- set up, instrument (after storage or transport) 73
- Shutter function test 57
- specification
 - halogen lamp 63

- installation 47
- instrument fuses 66
- spectra
 - CY3® dye 41
 - CY5® dye 42
 - FAM™ dye 41
 - JOE™ dye 41
 - NED™ dye 41
 - ROX™ dye 42
 - SYBR® Green I dsDNA Binding dye 41
 - TAMRA™ dye 41
 - TEXAS RED® dye 42
 - VIC® dye 41
- standard plate, description 9
- standards
 - EMC xxii
 - safety xxii
- storage, prepare instrument for inactivity 70
- SYBR® Green I dsDNA Binding dye 33, 41
- symbols, safety xiii
- system dye
 - CY3® dye 33, 41
 - CY5® dye 33, 42
 - FAM™ dye 33, 41, 46
 - JOE™ dye 33, 41
 - NED™ dye 33, 41
 - ROX™ dye 33, 42
 - SYBR® Green I dsDNA Binding dye 33, 41
 - TAMRA™ dye 33, 41
 - TEXAS RED® dye 33, 42
 - VIC® dye 33, 41

T

- TAMRA™ dye 33, 41
- TaqMan® Fast Universal PCR Master Mix, No AmpErase® UNG 46
- Technical Support, contacting x
- TEXAS RED® dye 33, 42
- Thermal Cycler function test 57
- training, information on x
- troubleshoot
 - background calibration 30
 - dye calibration 43
 - function tests 57
 - instrument fuses 66–67
 - instrument log 56
 - lamp replacement 63–65
 - lamp status, view 58
 - optical calibration 30
 - RNase P experiment 54
 - ROI calibration 17
 - sample block decontamination 59

U

- update
 - operating system 67

service packs 67
USB function test 57
user attention words, described vii

V

VIC[®] dye 33, 41

W

WARNING, description xii
waste disposal, guidelines xix
workstation safety xxi

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