

© Copyright 2006, 2010 Applied Biosystems. All rights reserved.

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

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:

The Applied Biosystems 7300, 7500 and 7500 Fast Real-Time PCR Systems are real-time thermal cyclers 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.

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

### TRADEMARKS:

AB (Design), Applied Biosystems, ABI PRISM, NucPrep, PrepMan, Primer Express, and VIC are registered trademarks and Applera, BloodPrep, FAM, MicroAmp, ROX, and TAMRA are trademarks of Applied Biosystems or its subsidiaries in the U.S. and/or other countries.

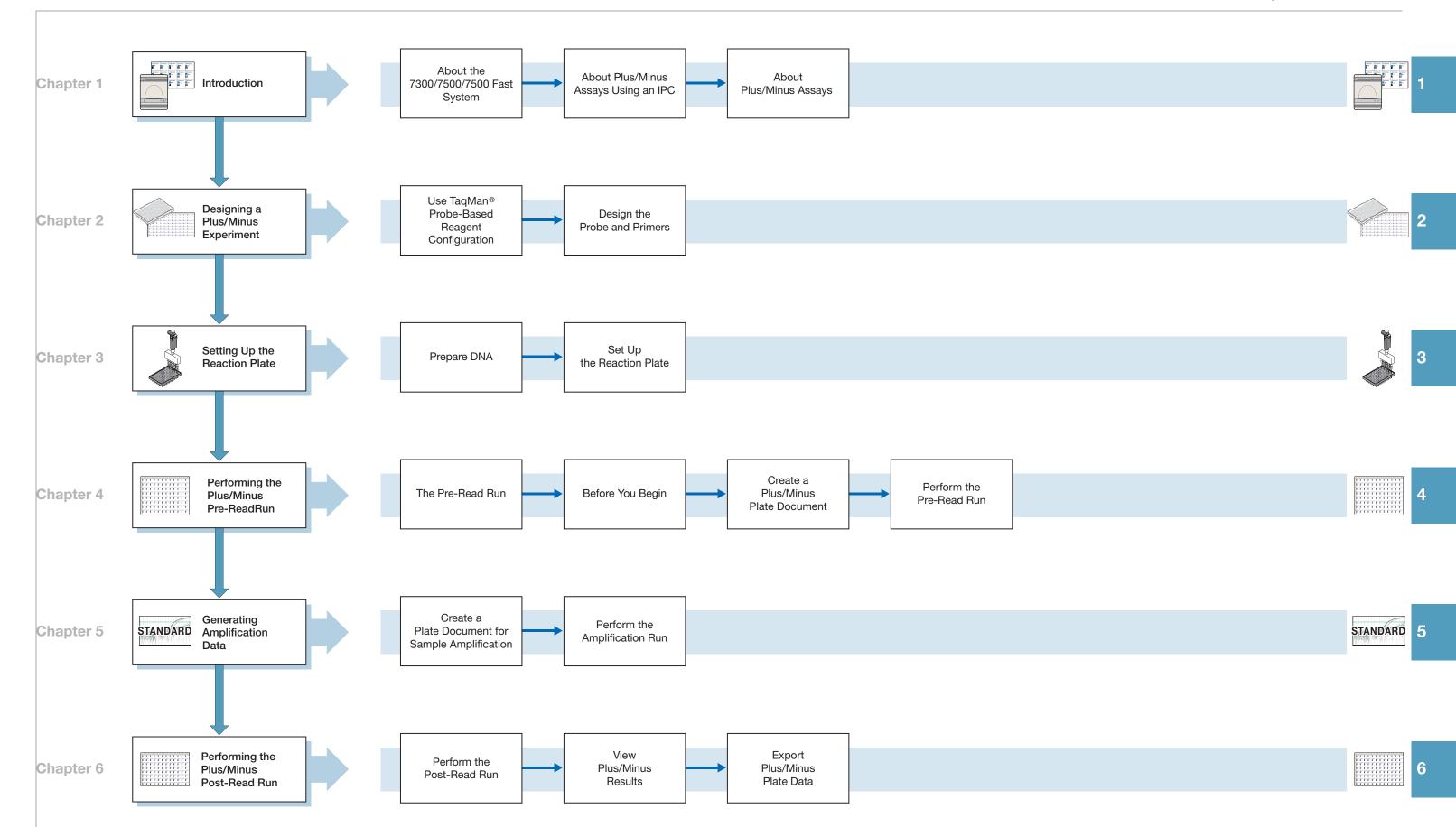
AmpErase, AmpliTaq Gold, and TaqMan are registered trademarks of Roche Molecular Systems, Inc.

SYBR Green is a registered trademark of Molecular Probes, Inc.

Excel, Microsoft, PowerPoint, and Windows are registered trademarks of Microsoft Corporation.

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

Part Number 4347821 Rev. E 06/2010



# Contents

	Plus/Minus Experiment Workflow	III
	Preface	vii
	How to Use This Guide	vii
	How to Obtain More Information	viii
	How to Obtain Support	ix
	Safety	хi
	Safety Alert Words	xi
	Good Laboratory Practices	xii
	General Chemical Warnings	
	General Biohazard Warnings	
	General Chemical Waste Hazard Warnings	
	Obtaining Modos	
Chapter 1	Introduction	1
	About the 7300/7500/7500 Fast System	2
	About Plus/Minus Assays Using an IPC	
	About Plus/Minus Assays	4
Chapter 2	Designing a Plus/Minus Experiment	7
	Using TaqMan® Probe-Based Reagent Configuration	8
	Designing the Probe and Primers	
Chapter 2	Satting Up the Depation Dieto	11
Chapter 3	Setting Up the Reaction Plate	
	Preparing DNA	
	Setting Op the neaction Flate	
Chapter 4	Performing the Plus/Minus Pre-Read Run	17
-	The Pre-Read Run	18
	Before You Begin	18
	Creating a Plus/Minus Plate Document	18

	Performing the Pre-Read Run	24
Chapter 5	Generating Amplification Data	27
	Creating a Plate Document for Sample Amplification	
Chapter 6	Performing the Plus/Minus Post-Read Run	35
	Performing the Post-Read Run	39
Appendix A	Creating Detectors	47
Appendix B	Viewing Amplification Data  Configuring Analysis Settings	
	Viewing the Amplification Data	
Appendix C	Example Plus/Minus Experiment	57
	References	65
	Index	67

## **Preface**

## How to Use This Guide

# Purpose of This Guide

This manual is written for principal investigators and laboratory staff who run plus/minus assays using the Applied Biosystems 7300/7500/7500 Fast Real-Time PCR System (7300/7500/7500 Fast system).

### Assumptions

This guide assumes that you have:

- Familiarity with Microsoft® Windows® XP operating system.
- Knowledge of general techniques for handling DNA samples and preparing them for PCR.
- A general understanding of hard drives and data storage, file transfers, and copying and pasting.

### **Text Conventions**

This guide uses the following conventions:

- Bold indicates user action. For example:
   Type 0, then press Enter for each of the remaining fields.
- *Italic* text indicates new or important words and is also used for emphasis. For example:

Before analyzing, *always* prepare fresh matrix.

 A right arrow bracket (>) separates successive commands you select from a dropdown or shortcut menu. For example:

Select File > Open.

# User Attention Words

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

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

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

**CAUTION** Indicates a potentially hazardous situation that, if not avoided, may result in minor or moderate injury. It may also be used to alert against unsafe practices.

**WARNING** Indicates a potentially hazardous situation that, if not avoided, could result in death or serious injury.

## **How to Obtain More Information**

# Related Documentation

For more information about using the 7300/7500/7500 Fast system, refer to:

- Applied Biosystems 7300/7500/7500 Fast Real-Time PCR System Online Help
- Applied Biosystems 7300/7500/7500 Fast Real-Time PCR System Absolute Quantitation Using Standard Curve Getting Started Guide (PN 4347825)
- Applied Biosystems 7300/7500/7500 Fast Real-Time PCR System Allelic Discrimination Getting Started Guide (PN 4347822)
- Applied Biosystems 7300/7500/7500 Fast Real-Time PCR System Relative Quantitation Using Comparative C<sub>T</sub> Getting Started Guide (PN 4347824)
- Applied Biosystems 7300/7500/7500 Fast Real-Time PCR System: User Guide for the 21 CFR Part 11 Module in SDS Software v1.4 (PN 4374432)
- Applied Biosystems 7300/7500/7500 Fast Real-Time PCR System Installation and Maintenance Guide (PN 4347828)
- Applied Biosystems 7300/7500/7500 Fast Real-Time PCR System Site Preparation Guide (PN 4347823)
- Applied Biosystems 7300/7500/7500 Fast Real-Time PCR System Performing Fast Gene Quantitation Quick Reference Card (PN 4362285)
- Applied Biosystems Real-Time PCR Systems Computer Setup Guide (PN 4365367)
- Applied Biosystems Real-Time PCR Systems Chemistry Guide (PN 4348358)

# Accessing the Online Help

Access the Online Help System by clicking ? in the tool bar of the SDS software window, or by selecting **Help > Contents and Index**.

### Send Us Your Comments

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

techpubs@appliedbiosystems.com

## **How to Obtain Support**

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

At the Support page, you can:

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

### Preface

How to Obtain Support

# Safety

## 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, 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*.

### **Examples**

The following examples show the use of safety alert words:

**IMPORTANT!** You must create a separate a Sample Entry Spreadsheet for each 96-well microtiter plate.

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

WARNING CHEMICAL HAZARD. Formamide. Exposure causes eye, skin, and respiratory tract irritation. It is a possible developmental and birth defect hazard. 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.

## **Good Laboratory Practices**

## PCR Good Laboratory Practices

PCR assays require special laboratory practices to avoid false positive amplifications (Kwok and Higuchi, 1989). The high throughput and repetition of these assays can lead to amplification of a single DNA molecule (Saiki et al., 1985; Mullis and Faloona, 1987).

- Wear a clean lab coat (not previously worn while handling amplified PCR products or used during sample preparation) and clean gloves when preparing samples for PCR amplification.
- Change gloves whenever you suspect that they are contaminated.
- Maintain separate areas, dedicated equipment, and supplies for:
- Sample preparation and PCR setup
- PCR amplification and post-PCR analysis
- Never bring amplified PCR products into the PCR setup area.
- Open and close all sample tubes and reaction plates carefully. Do not splash or spray PCR samples.
- Keep reactions and components sealed as much as possible.
- Use positive displacement pipettes or aerosol-resistant pipette tips.
- Clean lab benches and equipment periodically with freshly diluted 10% bleach solution.

## **Bibliography**

Kwok, S. and Higuchi, R. 1989. Avoiding false positives with PCR. *Nature* 339:237-238

Mullis, K.B. and Faloona, F.A. 1987. Specific synthesis of DNA *in vitro* via a polymerase-catalyzed chain reaction. *Methods Enzymol*. 155:335-350.

Saiki, R.K., Scharf, S., Faloona, F., *et al.* 1985. Enzymatic amplification of  $\beta$ - globin genomic sequences and restriction site analysis for diagnosis of sickle cell anemia. *Science* 230:1350-1354.

## **General Chemical Warnings**

## Chemical Hazard Warning

WARNING CHEMICAL HAZARD. Some of the chemicals used with Applied Biosystems instruments and protocols are potentially hazardous and can cause injury, illness, or death.

### Chemical Safety Guidelines

To minimize the hazards of chemicals:

- Read and understand the Material Safety Data Sheets (MSDS) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials.
- 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.

## General Warnings

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

## **General Biohazard Warnings**

## 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; <a href="http://bmbl.od.nih.gov">http://bmbl.od.nih.gov</a>)
- Occupational Safety and Health Standards, Bloodborne Pathogens (29 CFR§1910.1030; <a href="http://www.access.gpo.gov/nara/cfr/">http://www.access.gpo.gov/nara/cfr/</a> 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:

http://www.cdc.gov

## **General Chemical Waste Hazard Warnings**

## Chemical Waste Hazard

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

**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.

## **Obtaining MSDSs**

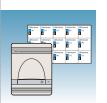
Chemical manufacturers supply current Material Safety Data Sheets (MSDSs) with shipments of hazardous chemicals to *new* customers. They also provide MSDSs with the first shipment of a hazardous chemical to a customer after an MSDS has been updated. MSDSs provide the safety information you need to store,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

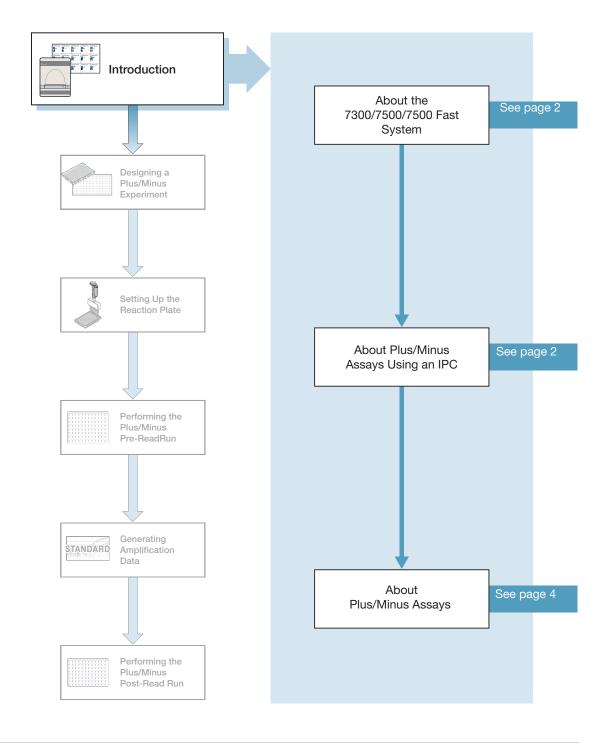
You can obtain from Applied Biosystems the MSDS for any chemical supplied by Applied Biosystems. This service is free and available 24 hours a day.

To obtain MSDSs:

- 1. Go to https://docs.appliedbiosystems.com/msdssearch.html
- 2. In the Search field, type in the chemical name, 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
- **4.** To have a copy of a document sent by fax or e-mail:
  - a. Select Fax or Email to the left of the document title in the Search Results page
  - **b.** Click **RETRIEVE DOCUMENTS** at the end of the document list.
  - c. After you enter the required information, click View/Deliver Selected Documents Now.



# Introduction



## About the 7300/7500/7500 Fast System

### Description

The Applied Biosystems Real Time PCR System (7300/7500/7500 Fast system) uses fluorescent-based PCR chemistries to provide:

- Quantitative detection of nucleic acid sequence using real-time analysis.
- Qualitative detection of nucleic acid sequence using end-point and dissociationcurve analysis.

### Plus/Minus Assay

The 7300/7500/7500 Fast system allows you to perform several assay types with plates in the 96-well format. This guide describes the plus/minus assay, which determines whether or not a specific target sequence is present in a sample.

**Note:** For information about the other assay types, refer to the *Real-Time PCR Systems Chemistry Guide* (PN 4348358) and the Online Help for the 7300/7500/7500 Fast system.

**Note:** Plus/Minus Assays may be run on a 7500 Fast system using standard reagents; Plus/Minus Assays are not supported using Fast reagents and protocols.

## About Plus/Minus Assays Using an IPC

### Definition

A plus/minus assay is an end-point assay that determines if a specific target sequence is present (plus) or not present (minus) in a sample. In an end-point assay, data are collected at the end of the PCR process.

### What Is An IPC?

An internal positive control (IPC) is used in plus/minus assays to monitor the PCR process and to ensure that a negative result is not due to failed PCR (see TaqMan® Exogenous Internal Positive Control Reagents, PN 4308323). The IPC consists of a template, a primer set, and a dye-labeled (VIC®) probe added to each well of a reaction plate. The IPC is part of the reaction mix. See "Preparing the PCR Reaction Mix" on page 14.

Plus/Minus assays with an IPC use fluorogenic 5' nuclease chemistry (TaqMan probebased chemistry). During amplification, the sample target and the IPC target generate reporter fluorescence signals, so that positive or negative calls may be made on unknown samples.

**Note:** The SYBR® Green I dye chemistry is not supported for plus/minus assays using an IPC.

Notes			



# Terms Used in Plus/Minus Analysis

Term	Definition
Internal positive control (IPC)	A second TaqMan <sup>®</sup> probe and primer set added to the reaction plate to monitor the PCR process and to ensure that a negative result is not due to failed PCR in the sample
No amplification control (NAC)	Wells that contain no target template and blocked IPC (the IPC target template has been blocked by a blocking agent)
No template control (NTC)	A sample that contains no target template
Nucleic acid target	Nucleotide sequence that you want to identify as present or absent
Unknown sample (U)	The sample for which you want to determine the presence or absence of a specific target

Notes\_\_\_\_

## **About Plus/Minus Assays**

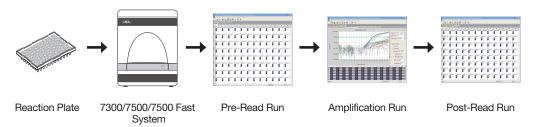
## Plus/Minus Experiment Workflow

This document uses the term "plus/minus assay" to refer to the entire process of analyzing samples of extracted DNA from data collected at the end of the PCR process.

Design the experiment and isolate the DNA, then conduct a plus/minus assay by performing:

- **A pre-read run** on a plus/minus plate document to determine the baseline fluorescence associated with primers and probes before amplification.
- An amplification run using an AQ plate document to generate real-time PCR data, which can be used to analyze and troubleshoot the PCR data for the plus/minus assay, if needed.
- A post-read run using the original plus/minus plate document, which automatically subtracts the baseline fluorescence determined during the pre-read run to calculate the result.

The following figure illustrates the complete process.



## Required User-Supplied Materials

lå	Sauras
Item	Source
DNA isolation and purification chemistry systems:	
ABI PRISM® 6100 Nucleic Acid PrepStation	Applied Biosystems (PN 6100-01)
<ul> <li>BloodPrep<sup>™</sup> Chemistry (genomic DNA from fresh or frozen blood or cells)</li> </ul>	Applied Biosystems (PN 4346860)
<ul> <li>NucPrep<sup>®</sup> Chemistry (DNA from animal and plant tissue)</li> </ul>	Applied Biosystems (PN 4340274)
PrepMan® Ultra Sample Preparation Reagent with Protocol	Applied Biosystems (PN 4322547)
Labeled primers and probes source:	
Primer Express® Software (custom-	• PN 4363991 (1-user license)
designed primers and probes)	• PN 4363993 (5-user license)
MicroAmp <sup>™</sup> Optical 96-Well Reaction Plates with Barcode	Applied Biosystems (PN 4306737)
MicroAmp <sup>™</sup> Optical Adhesive Film	Applied Biosystems (PN 4311971)
6700 Reagent Tubes, 10-mL	Applied Biosystems (PN 4305932)

Notes		

Item	Source
TaqMan <sup>®</sup> Exogenous Internal Positive Control Reagents	Applied Biosystems (PN 4308323)
TaqMan <sup>®</sup> Universal PCR Master Mix	Applied Biosystems (PN 4304437)
Centrifuge with adapter for 96-well plates	Major Laboratory Supplier (MLS)
Gloves	MLS
Microcentrifuge	MLS
Microcentrifuge tubes, sterile 1.5-mL	MLS
Nuclease-free water	MLS
Pipette tips, with filter plugs	MLS
Pipettors, positive-displacement	MLS
Tris-EDTA (TE) Buffer, pH 8.0	MLS
Vortexer	MLS

## **Example Plus/Minus Experiment**

To better illustrate how to design, perform, and analyze plus/minus experiments, this document provides an example experiment. The example experiment represents a typical plus/minus experiment that you can use as a quick-start procedure to familiarize yourself with the plus/minus workflow. Details about the plus/minus workflow are described in the subsequent chapters of this guide. Example Experiment boxes appear in subsequent chapters to illustrate workflow details. Refer to Appendix C, Example Plus/Minus Experiment on page 57 for more information. To view the example experiment data file in the SDS software:

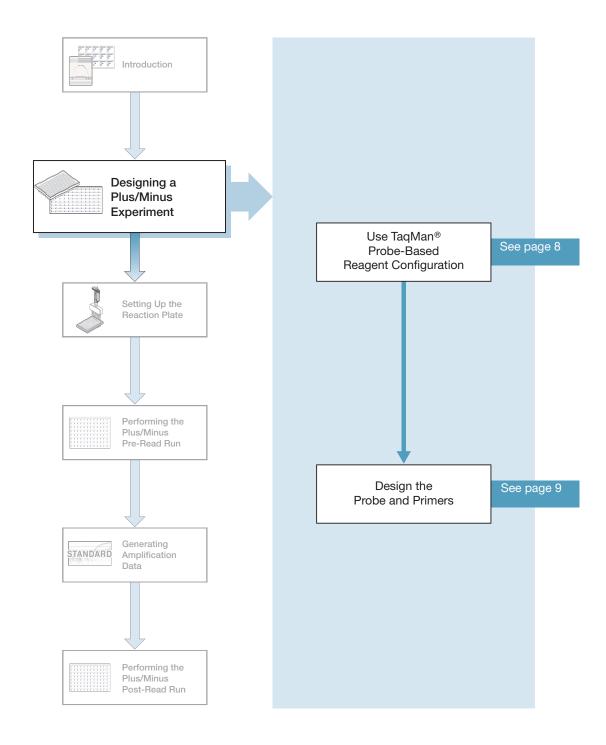
- 1. Select File > Open.
- 2. Navigate to Applied Biosystems\SDS Documents\Example Data Files\ EXAMPLE PM.sds, then click Open.



Notes	
6	Plus/Minus Assay Getting Started Guide for 7300/7500/7500 Fast Systems



# Designing a Plus/Minus Experiment



# Using TaqMan® Probe-Based Reagent Configuration

## About the Chemistry

Plus/minus assays with an IPC use the fluorogenic 5' nuclease chemistry (TaqMan<sup>®</sup> probe-based chemistry) to detect a specific PCR product as it accumulates during PCR cycles. For more information about the TaqMan probe-based chemistry, refer to the *Real-Time PCR Systems Chemistry Guide* (PN 4348358).

## PCR and Detection of cDNA a. Assay Components Reverse primer Forward primer **LEGEND** Random Primer b. Denatured Template and Annealing of Assay Components Reverse Transcriptase FAM™ dye Forward prime Quencher Minor Groove c. Signal Generation AmpliTaq Gold® DNA Polymerase Probe Primer Template Extended Primer

Notes			



# Chemistry Kits for Plus/Minus Assay

The following reagents are available from Applied Biosystems for designing and running plus/minus assays.

Kit	Part Number
TaqMan® Exogenous Internal Positive Control Reagents	4308323
TaqMan® Universal PCR Master Mix	4304437

**Note:** The IPC DNA, primers, and probe supplied in these reagents can be used with all sample target systems. Refer to the *TaqMan*<sup>®</sup> *Universal PCR Master Mix Protocol* (PN 4304449) for instructions on optimizing amplification of your target.

## **Designing the Probe and Primers**

Design a probe and primer set for your target sequence. Applied Biosystems provides the Primer Express® software for this purpose. For more information about using this software, refer to the *Primer Express Software v3.0 Getting Started Guide* (PN 4362460).

### **Example Experiment**

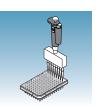
To conduct the example experiment, DNA was extracted from 84 batches of hamburger meat and tested for the presence of *E. coli* using the plus/minus assay on the 7300 Real Time PCR System. Six no IPC/no target template controls, six IPC/no target template controls, and 84 unknown samples were run.

The TaqMan<sup>®</sup> Exogenous Internal Positive Control Reagents supplied one 1-mL tube of 10X Exo IPC Mix. This mix contained the IPC primers and VIC<sup>®</sup>-labeled probe. The primers/probe set for *E. coli* was designed using Applied Biosystems Primer Express software, and contained a FAM<sup>™</sup> dye-labeled probe with a TAMRA<sup>™</sup> dye-labeled quencher.

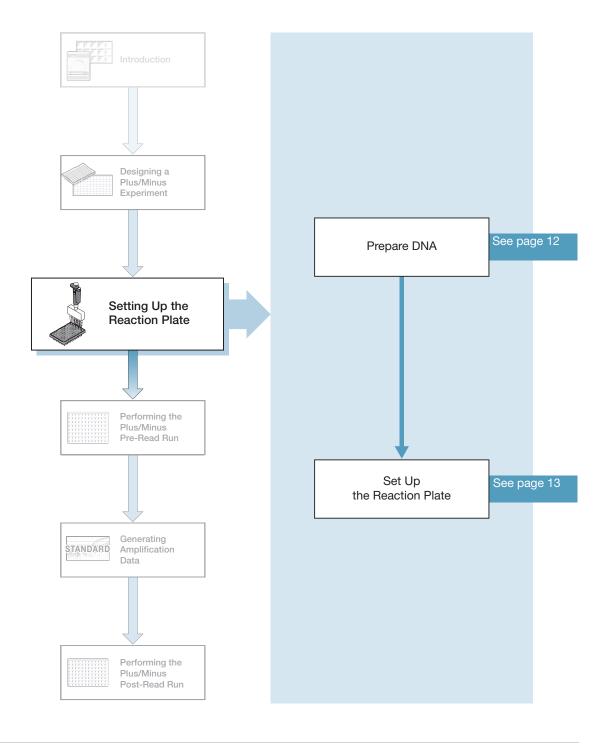
Notes		



Notes	
10	Plus/Minus Assay Getting Started Guide for 7300/7500/7500 Fast Systems



# Setting Up the Reaction Plate



## **Preparing DNA**

## Systems and Chemistries for DNA Isolation

Applied Biosystems supplies several instrument systems and chemistries for isolating DNA from a variety of starting materials, such as blood, tissue, cell cultures, plant material, and food.

System	Part Number
BloodPrep <sup>™</sup> Chemistry	4346860
NucPrep <sup>®</sup> Chemistry	4340274
PrepMan® Ultra Sample Preparation Reagent with Protocol	4322547
ABI PRISM® 6100 Nucleic Acid PrepStation	6100-01

For more information, refer to:

- DNA Isolation from Fresh and Frozen Blood, Tissue Culture Cells, and Buccal Swabs Protocol (PN 4343586)
- NucPrep® Chemistry Isolation of Genomic DNA from Animal and Plant Tissue Protocol (PN 4333959)

### **Quality of DNA**

Ensure that the DNA you use for a plus/minus experiment:

- Has a A<sub>260/280</sub> ratio of greater than 1.7.
- Is extracted from the raw material you are testing using an optimized protocol.
- Does not contain PCR inhibitors.
- Is intact as visualized by gel electrophoresis.
- Has not been heated above 60 °C, which can cause degradation.

### **Example Experiment**

The meat samples were frozen with liquid nitrogen and ground to a fine powder with a pre-chilled mortar and pestle. DNA was extracted using the PrepMan<sup>®</sup> Ultra Sample Preparation Reagent with Protocol (PN 4322547) to obtain a final concentration of 10  $ng/\mu L$  of DNA for each sample.

Notes		



## **Setting Up the Reaction Plate**

This section describes how to set up a 96-well plate for a plus/minus run with samples and reaction mix. The reagents, volumes, and final concentrations in "Preparing the PCR Reaction Mix" on page 14 were taken from the *TaqMan*® *Exogenous Internal Positive Control Reagents Protocol* (PN 4308323).

## **Example Experiment** Extracted DNA samples were pipetted onto a 96-well plate along with negative and positive controls. Wells A1-A6 contained blocked IPC and no target template; wells A7-A12 contained IPC template (IPC+), but no target template; and wells B1-H12 contained both IPC and target template. Blocked IPC and no target 5 10 11 12 NAC NTC IPC and no target В C D E Unknown plus IPC IPC plus target G

Notes			

## Preparing the PCR Reaction Mix

Make a volume of reaction mix sufficient to provide 45  $\mu$ L (18  $\mu$ L for the Fast system) for each well you use on the plate.

CAUTION CHEMICAL HAZARD. TaqMan Universal PCR Master Mix (2X) No AmpErase UNG may cause eye and skin irritation. Exposure may cause discomfort if swallowed or inhaled. Read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Item	Volume for one Reaction (Standard)	Volume for one Reaction (Fast)	Final Concentration
TaqMan <sup>®</sup> Universal PCR Master Mix (2X)	25.0	10.0	1X
10X Exo IPC Mix (IPC kit)	5.0	2.0	50 to 900 nM
50X Exo IPC DNA (IPC kit)	1.0	0.4	50 to 900 nM
Target primers, probe, and deionized water	14.0	5.6	50 to 250 nM
Total	45.0	18.0	_

Notes\_\_\_\_

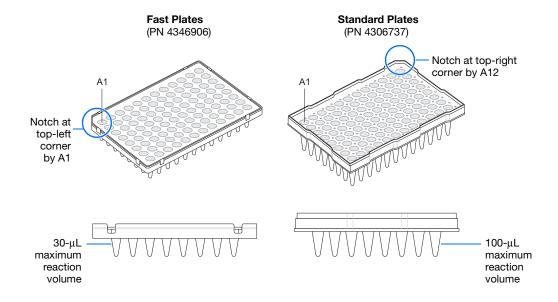


## Preparing the Plate

### Standard vs. Fast Plates

**IMPORTANT!** Ensure you use the standard Optical 96-Well Plate on the 7500 Real-Time PCR system. Fast Optical 96-Well Plates will *not* fit into the standard block correctly and using them will result in loss of data.

**IMPORTANT!** Ensure you use the Fast Optical 96-Well Plate on the 7500 Fast Real-Time PCR system. Standard plates will not function properly and they may be crushed when using the 96-Well Fast Block.



## To prepare the plate

- 1. Into each well, pipette 45  $\mu$ L (18  $\mu$ L for the Fast system) of the reaction mix.
- 2. Pipette 5 μL (2 μL for the Fast system) of sample (NAC, NTC, or unknowns) into each well of a 96-well plate:

	1	2	3	4	5	6	7	8	9	10	11	12
Α			N	AC					N	C		
В												
С												
D												
Ε						Sam	ples					
F					nkn	own	plu	SIP				
G												
Н												

**Note:** The final reaction volume in each well should be  $50 \mu L$  ( $20 \mu L$  for the Fast system).

**3.** Keep the reactions on ice until the plate is loaded into the 7300/7500/7500 Fast instrument.

If preparing	Then add (Standard)	Then add (Fast)
NAC <sup>‡</sup>	5 μL of 10X Exo IPC Block (IPC kit)	2 μL of 10X Exo IPC Block (IPC kit)
NTC§	5 μL of 1X TE or H <sub>2</sub> O	2 μL of 1X TE or H <sub>2</sub> O

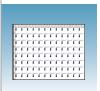
- No Amplification Control well contains no target template and blocked IPC.
- § No Template Control well contains no target template, only IPC.

### **Example Experiment**

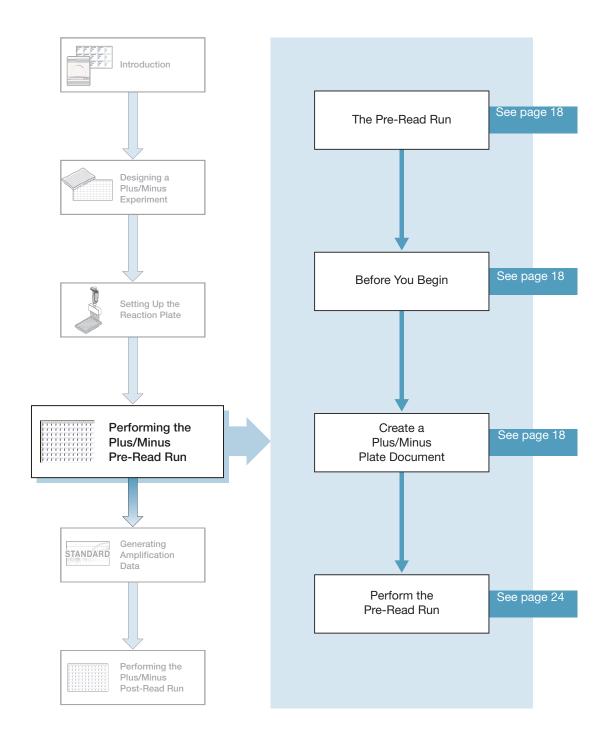
 $45~\mu\text{L}$  of reaction mix was pipetted into each well of a 96-well plate, and  $5~\mu\text{L}$  of the following was added as specified in the table below.

Wells	To prepare	Add to each well (Standard)	Add to each well (Fast)
A1 to A6	NAC‡	5 μL of 10X Exo IPC Block (IPC kit)	2 μL of 10X Exo IPC Block (IPC kit)
A7 to A12	NTC§	5 μL of 1X TE or H <sub>2</sub> O	2 μL of 1X TE or H <sub>2</sub> O
B1 to H12	U#	5 μL of sample DNA	2 μL of sample DNA

- § No Template Control well contains no target template, only IPC.
- # Unknown well contains both target template and IPC.



# Performing the Plus/Minus Pre-Read Run



## The Pre-Read Run

A pre-read run records the background fluorescence of each well of the plus/minus plate before PCR. During the post-read run, the pre-read fluorescence is subtracted from the post-read fluorescence to account for pre-amplification background fluorescence. Subtracting pre-read from post-read fluorescence ensures accurate results.

## **Before You Begin**

Check that background and pure-dye runs have been performed regularly to ensure optimal performance of the 7300/7500/7500 Fast system. For more information about calibrating the system, see the *Applied Biosystems* 7300/7500/7500 Fast Real-Time PCR System Online Help.

## Creating a Plus/Minus Plate Document

A plus/minus plate document is an SDS software document that stores data collected from a plus/minus run for a single 96-well plate. Plus/Minus plate documents also store other information about the run, including sample names and detectors.

### Plate Document Parameters

When you create a plus/minus plate document with an IPC, you define specific parameters for each plus/minus reaction plate:

- **Detectors** A virtual representation in the SDS software of a TaqMan<sup>®</sup> probe and primer set and an associated fluorescent dye that detects a single target nucleic acid sequence. Appendix A on page 47 explains how to create detectors.
- Task A setting that you apply to each well of a plate document that determines the way the SDS software uses the data collected from the well during analysis.

**Note:** Applied Biosystems recommends that you run six replicates of each control (NAC and NTC) to accurately define plus/minus thresholds and to obtain plus/minus calls with a 99.7% confidence level.

Notes		



## **Detector Tasks**

You assign a task to each detector in each well of a plate document.

For plus/minus plate documents, there are four types of tasks:

Task	Symbol	Apply to
Unknown	U	All detectors of wells that contain target sequence.
IPC	1	All detectors of wells that contain IPC.
IPC+	I <del>I</del>	All detectors of control wells that contain IPC with a blocking agent but no target template.
NTC	N	All detectors of negative control wells that contain PCR reagents, but no target template and no IPC.

# Creating a New Plus/Minus Plate Document

You can enter sample information into a new plate document, copy or import sample information from existing plate documents, or use a template document to set up new plate documents.

This section describes setting up new plate documents. Refer to the Online Help for information about copying or importing sample information or using template documents.

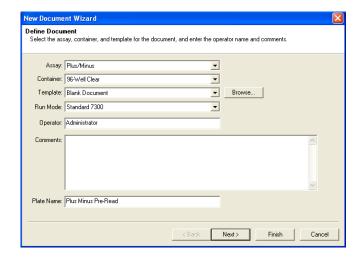
**Note:** The following procedure is illustrated using the example experiment data file (see page 5).

### To create a new plus/minus plate document:

- 1. Select Start > All Programs > Applied Biosystems > 7300/7500/7500 Fast System > 7300/7500/7500 Fast System Software () to start the 7300/7500/7500 Fast SDS software.
- 2. In the Quick Startup document dialog box, select Create New Document.



- 3. In the New Document Wizard, click the assay drop-down list, then select Plus/Minus assay. Accept the default settings for the Container and Template fields (96-Well Clear and Blank Document).
- **4.** In the Plate Name field, type **Plus Minus Pre-Read**.
- 5. Click Next>.





- **6.** Select the detectors to add to the plate document.
  - **a.** Click to select a detector. (Ctrl-click to select multiple detectors.)

**Note:** If no detectors are listed, click **New Detector** to open the New Detector dialog box. For more information about creating new detectors, refer to Appendix A, "Creating Detectors," on page 47.

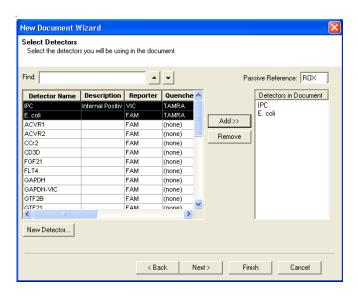
**b.** Click **Add>>**. The detector(s) are added to the Detectors in Document list box.

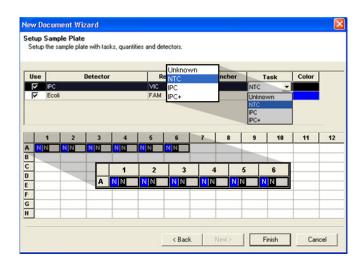
**IMPORTANT!** Ensure that the reporter dye for the target is different from the reporter dye for the IPC, which is VIC® dye.

**Note:** The TaqMan<sup>®</sup> Exogenous Internal Positive Control Reagents (PN 4308323) use an IPC VIC dye-labeled probe with TAMRA<sup>™</sup> dyelabeled quencher.

**Note:** To remove a detector in the Detectors in Document window, select the detector, then click **Remove**.

- c. Click Next.
- **7.** Select six wells on the plate document for the no amplification controls (blocked IPC and no target template).
  - **a.** Select wells A1 to A6.
  - b. Select the target detector by checking the Use box next to it.
  - c. Select Task > NTC. An N appears in the well.
  - d. Select the IPC detector by checking the Use box next to it.
  - e. Select Task > NTC. A second N appears in the well.







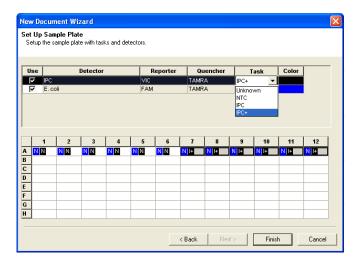
- **8.** Select six more wells on the plate document for the no template controls (IPC but no target template).
  - a. Select wells A7 to A12.
  - b. Select the target detector by checking the Use box next to it.
  - c. Select Task > NTC. An N appears in the well
  - d. Select the IPC detector by checking the Use box next to it.
  - e. Select **Task > IPC**<sup>+</sup>. An I+ appears in the well next to the N.
- **9.** Select all remaining wells for the unknown samples (IPC and target template).
  - **a.** Select wells B1 to H12 by click-dragging across all empty wells.
  - **b.** Select the target detector by checking the Use box next to it.
  - **c.** Select **Task > Unknown**. A U appears in the boxes.
  - d. Select the IPC detector by checking the Use box next to it.
  - e. Select **Task > IPC**. An I appears next to the U.

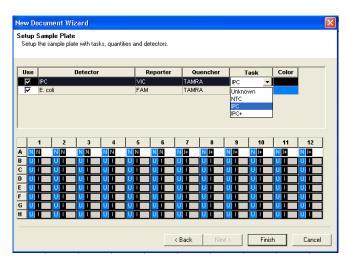
### 10. Click Finish.

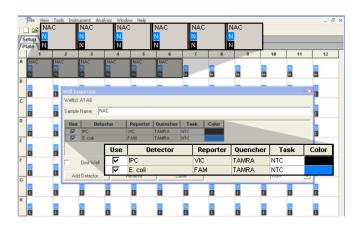
- **11.** Enter sample names for each well.
  - a. Click or select View > Well Inspector from the menu.

**Note:** To enter sample names without using the Well Inspector, click-drag to select wells, then type the sample name.

- **b.** Click-drag to select replicate wells.
- **c.** Type the sample name.







- d. Accept the default setting (ROX<sup>™</sup>) for the Passive Reference, ROX dye.
- **e.** Repeat steps b through c until you name all wells.
- f. Click I to close the Well Inspector.

**IMPORTANT!** If your experiment does not use all the wells on a plate, do not omit the wells from use at this point. You can omit unused wells after the run is completed. For more information about omitting wells, refer to Online Help.

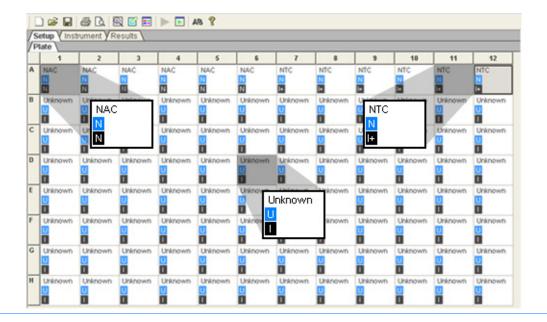
**Note:** You can change the sample setup information (sample name, detector, task) after a run is complete.

**g.** Verify the information on each well in the Setup tab.



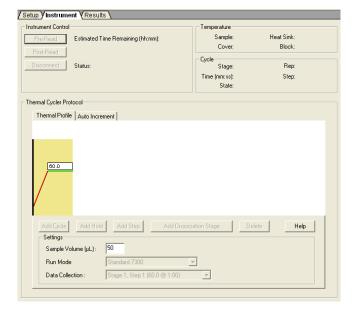
### **Example Experiment**

The pre-read plate document we created with controls and samples is shown in the picture below. We selected two detector tasks for each well, one for the target and one for the IPC. There are three possible combinations shown below: the NN box is the NAC (no target template, plus blocked IPC), the NI<sup>+</sup> box seen below is the NTC (no target template plus IPC), and the UI box is the unknown sample plus IPC.



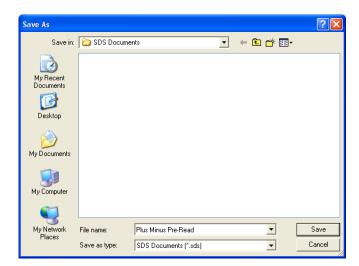
# Performing the Pre-Read Run

- 1. Select the **Instrument** tab.
- **2.** Accept the default value for sample volume.



3. Select File > Save, type Plus Minus Pre-Read for the plus/minus plate document, then click Save.

(Optional) If you want to use this plate document again, save it as a template document. Select File > Save As. In the Save in drop-down list, navigate to Applied Biosystems\7300\7500\7500 Fast System\Templates. Type a File name, then select Save as type: (\*.sdt).

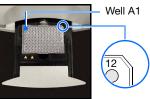




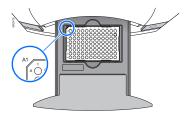
**4.** Load the reaction plate into the instrument.

**Note:** The A1 position is in the top-left side of the instrument tray.





7300/7500 system: Notched top-right corner for standard plates



7500 Fast system: Notched top-left corner for Fast plates

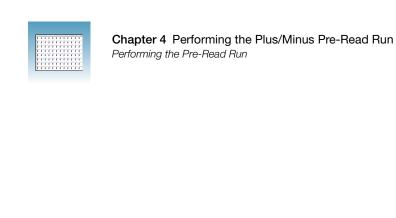
### 5. Click Pre-Read.

During the pre-read run, the instrument collects one fluorescence scan per well.

As the instrument performs the run, it displays status information in the Instrument tab. After the run is finished, the status values and the buttons are greyed out, and a message indicates whether or not the run is successful.

### 6. Select File > Close.

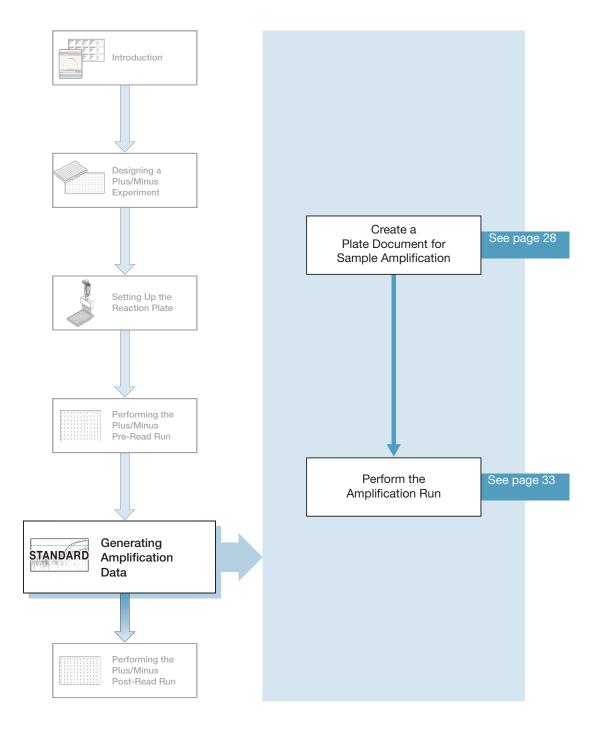




Notes\_\_\_\_



# **Generating Amplification Data**



## Creating a Plate Document for Sample Amplification

### Benefits of Real-Time Amplification

Because the plus/minus assay is an end-point assay, you can amplify the target sequences offline using any thermal cycler. However, using the 7300/7500/7500 Fast system to amplify the target sequences provides real-time PCR data. After the plus/minus samples are analyzed, you can study the amplification plots if you observe questionable calls or observe no data for a well.

### Using AQ Plate Documents for Amplification

You create and use Absolute Quantitation (AQ) plate documents to store real-time data for plus/minus assays. Because the AQ plate document is used only to amplify target sequences (not to quantify the PCR data), you do not need a standard curve for the AQ plate.

### **Detector Tasks**

For AQ plate documents, there are three types of tasks:

Task	Symbol	Apply to
Unknown	U	Detectors of wells that contain target sequences.
Standard	S	Detectors of wells that contain samples of known quantities.
NTC	N	Detectors of negative control wells that contain no template.

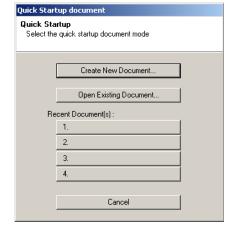
The task label "unknown" is used for both IPC and the target samples.

Notes		



### To create a new AQ plate document:

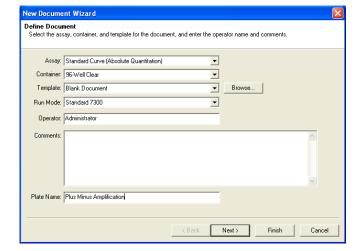
- 1. Select Start > All Programs > Applied Biosystems > 7300/7500/7500 Fast System > 7300/7500/7500 Fast System Software () to start the 7300/7500/7500 Fast SDS instrument software.
- 2. In the Quick Startup document dialog box, select Create New Document.



3. In New Document Wizard Assay drop-down list, select Standard Curve (Absolute Quantitation). Accept the default settings for the Container and Template (96-Well Clear and Blank Document).

**Note:** A standard curve is not necessary for a non-quantitation amplification run.

- **4.** In the Plate Name field, type **Plus Minus Amplification**.
- **5.** Click **Next>**. A list of detectors is displayed in the New Document Wizard.





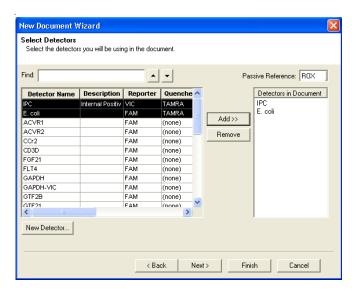
- **6.** Select the detectors to add to the plate document.
  - a. Click to select a detector. (Ctrl-click to select multiple detectors, for example IPC and E. coli.)
  - **b.** Click **Add>>.** The detector(s) are added to the Detectors in Document list box.

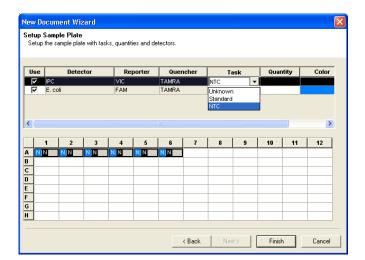
**IMPORTANT!** Ensure that the reporter dye for the target detector is different from the reporter dye for the IPC detector, which is VIC® dye.

Note: The TaqMan<sup>®</sup> Exogenous Internal Positive Control Reagents kit uses an IPC VIC dye-labeled probe with TAMRA<sup>™</sup> dye-labeled quencher.

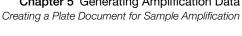
**Note:** To remove a detector in the Detectors in Document window, select the detector, then click **Remove**.

- c. Click Next.
- **7.** Select six wells on the plate document for the no amplification control (blocked IPC and no target template).
  - a. Select wells A1-A6.
  - b. Select the target detector by checking the Use box next to it.
  - c. Select Task > NTC. An N appears in the well.
  - d. Select the IPC detector by checking the Use box next to it.
  - e. Select Task > NTC. A second N appears in the well.





# Chapter 5 Generating Amplification Data



- **8.** Select six more wells on the plate document for the no template control (IPC but no target template).
  - a. Select wells A7-A12.
  - **b.** Select the target detector by checking the Use box next to it.
  - c. Select Task > NTC. An N appears in the well
  - d. Select the IPC detector by checking the Use box next to it.
  - e. Select Task > Unknown. A U appears in the well next to the N.
- **9.** Select all remaining wells for the unknown samples (IPC template and target template).
  - a. Select wells B1-H12 by click-dragging across all empty wells.
  - **b.** Select your **Target** detector by checking the Use box next to it.
  - c. Select Task > Unknown. A U appears in the boxes.
  - d. Select the IPC detector by checking the Use box next to it.
  - e. Select Task > Unknown. Another U appears next to the target U.

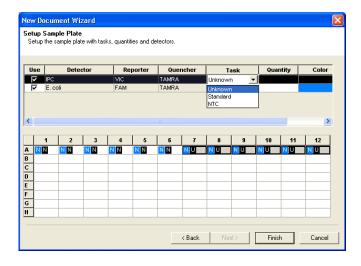
### 10. Click Finish.

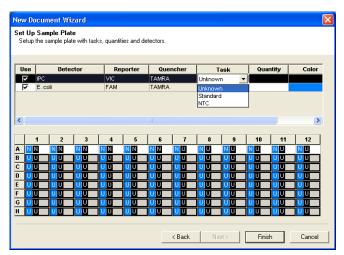
The 7300/7500/7500 Fast SDS software creates the plate document.

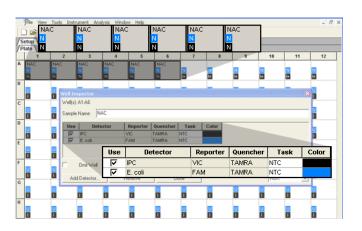
- **11.** Enter sample names for each well.
  - a. Double-click one well to open the Well Inspector, or select **View > Well Inspector**.

**Note:** To enter sample names without using the Well Inspector, click-drag to select wells, then type the sample name.

**b.** Click-drag to select all replicate wells for that sample.







### Chapter 5 Generating Amplification Data Creating a Plate Document for Sample Amplification



- **c.** Type the sample name in the Well Inspector. The information appears in the selected well(s).
- d. Accept the default setting (ROX<sup>™</sup> dye) for the Passive Reference, ROX dye. Optionally, you can change the detector task and Passive Reference dye.
- **e.** Repeat steps b through d until all wells have names.
- f. Click **▼** to close the Well Inspector.

**IMPORTANT!** If your experiment does not use all the wells on a plate, do not omit the wells from use at this point. You can omit unused wells after the run. For information about omitting unused wells, refer to the Online Help.

**g.** Verify the information about each well in the Setup tab.

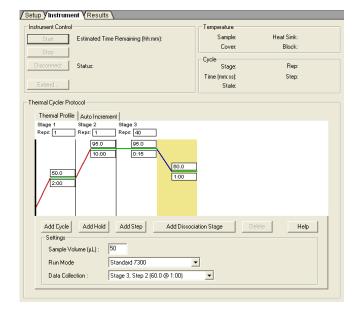
Notes\_\_\_\_\_



# Performing the Amplification Run

### 1. Select the **Instrument** tab.

By default, the standard PCR conditions for the PCR step are displayed.



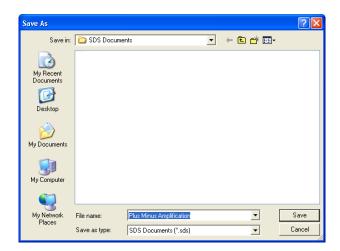
Times and Temperatures				
Initial Steps PCR (Each of 40 cycles)				
AmpErase® UNG Activation	AmpliTaq Gold <sup>®</sup> DNA Polymerase Activation	Melt	Anneal/Extend	
HOLD	HOLD	CYCLE	CYCLE	
2 min @ 50 °C	10 min @ 95 °C	15 sec @ 95 °C	1 min @ 60 °C	

### **2.** Verify the:

- Sample volume  $50 \mu L$  (20  $\mu L$  for the Fast system).
- Run Mode (for example, Standard 7300).



**3.** Select **File** > **Save**, then click **Save** to retain the name you assigned when you created the plate document.



**4.** Load the reaction plate into the instrument.

**Note:** The A1 position is in the top-left side of the instrument tray.

### 5. Click Start.

As the instrument performs the PCR run, it displays real-time status information in the Instrument tab and records the fluorescence resulting from cleavage of TaqMan® probes in the presence of the target sequences.

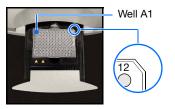
After the run is finished, the status values and the buttons are grayed-out, the Analysis button is enabled (), and a message indicates whether or not the run is successful.

All data generated during the run are saved to the plate document that you specified in step 3, and this data can be analyzed later for troubleshooting purposes.

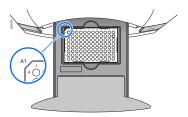
6. To view real-time PCR after the run is finished, click the Analysis button ▶, select the **Results** tab, select the **Amplification Plot** tab, then select all wells in the upper left box (next to A1).

**Note:** You can change the sample setup information (sample name, detector, task) after a run is complete.

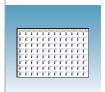




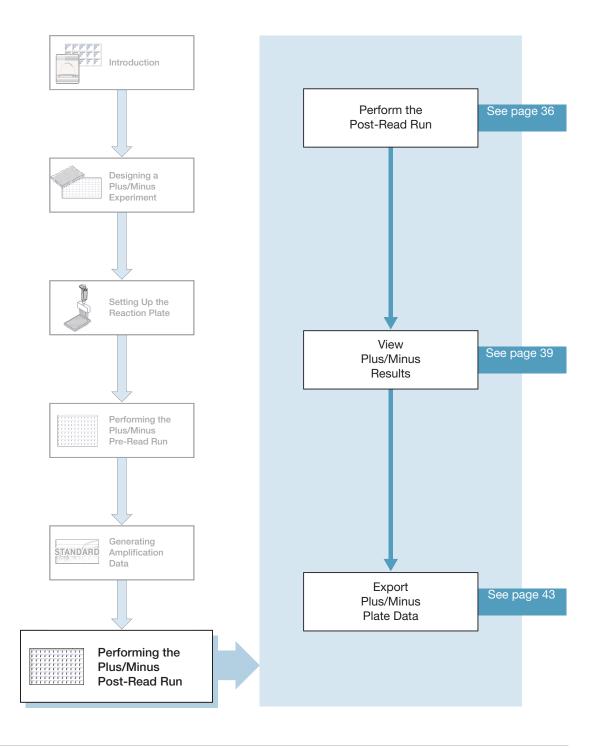
7300/7500 system: Notched top-right corner for standard plates



7500 Fast system: Notched top-left corner for Fast plates



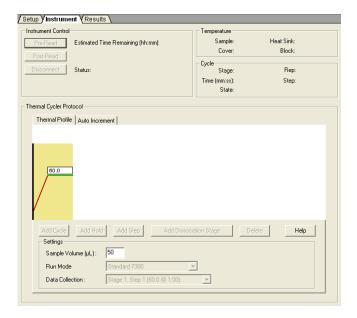
# Performing the Plus/Minus Post-Read Run



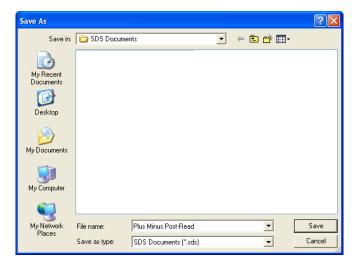
# Performing the Post-Read Run

Open the pre-read plate document.

1. Select the **Instrument** tab.



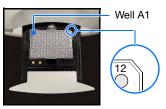
- **2.** Accept the default value for sample volume.
- 3. Select File > Save As, type the name Plus Minus Post-Read for the plus/minus plate document, then click Save.



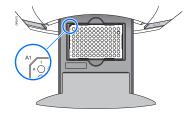
**4.** Load the reaction plate into the instrument.

**Note:** The A1 position is in the top-left side of the instrument tray.





7300/7500 system: Notched top-right corner for standard plates



7500 Fast system: Notched top-left corner for Fast plates

### 5. Click Post-Read.

After the run is finished, the status values and buttons are grayed-out, and a message indicates whether or not the run is successful.

**6.** Click the green analysis button (▶) to start analysis.

All data generated during the run are saved to the plus/minus plate document that you specified in step 3.

Notes

6





### **Example Experiment**

In the example plus/minus experiment using an IPC, the pre-read run was subtracted from the post-read run to account for background fluorescence. Post-read results for the presence of *E. coli* are displayed in the Results > Plate tab. For an explanation of results see "Viewing Plus/Minus Results" on page 39.

NAC	NAC N	NAC N	NAC N	NAC N	NAC N	NTC N	NTC N	NTC N	NTC N	NTC N	NTC N
	N	N	N	N	N	l+	l+	I <del>+</del>	I <del>+</del>	I <del>.</del>	<b>E</b>
Inknown (+)	Unknown U (+)	Unknown U (?)	Unknown U (-)	Unknown U (+)	Unknown U (-)						
nknown (+)	Unknown U (-)	Unknown U (+)	Unknown U (+)	Unknown U (+)	Unknown U (+)						
Inknown (+)	Unknown	Unknown U (+)	Unknown	Unknown U (+)	Unknown U (+)	Unknown U (+)	Unknown U (-)				
Inknown (+)	Unknown	Unknown U (+)	Unknown	Unknown U (+)	Unknown U (+)	Unknown U (+)	Unknown U (+)	Unknown U (-)	Unknown U (-)	Unknown	Unknown U (+)
nknown (+)	Unknown	Unknown U (+)	Unknown	Unknown U (+)	Unknown	Unknown	Unknown U (-)				
nknown (+)	Unknown	Unknown	Unknown U (+)	Unknown	Unknown	Unknown	Unknown	Unknown	Unknown	Unknown	Unknown U (+)
Inknown (+)	Unknown	Unknown U (-)	Unknown	Unknown	Unknown	Unknown	Unknown	Unknown	Unknown	Unknown	Unknown

Notes			

# **Viewing Plus/Minus Results**

### Results

After completing the plus/minus post-read run, the SDS software compares the relationship between the spectral changes in the unknown samples and the control reactions defined previously, NAC and NTC. An IPC threshold is calculated from the NAC control reactions, and the target threshold is calculated from the NTC control reactions. The target threshold is used to determine amplification of the unknown sample signal for each well. The IPC threshold is used to determine amplification of the IPC signal in each unknown sample well.

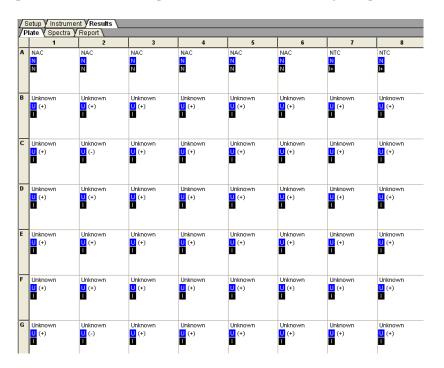
### Calling Unknowns Using IPC

To call the unknown samples, the SDS software compares the normalized reporter signal of each unknown sample to the target threshold and IPC sample to IPC threshold. The results are determined as follows:

- If the unknown sample signal is above the target threshold, then the call is positive (+).
- If the unknown sample signal is below the target threshold, the SDS software compares the IPC sample signal to the IPC threshold as follows:
  - If IPC sample signal is above the IPC threshold, then the call is negative (-).
  - If IPC sample signal is below the IPC threshold, then the call is undetermined (?).

### The Plate Tab

When the post-read is complete, select the Plate tab on the Results page to display the plus/minus calls for the presence or absence of the target sequence for each well.



Notes		

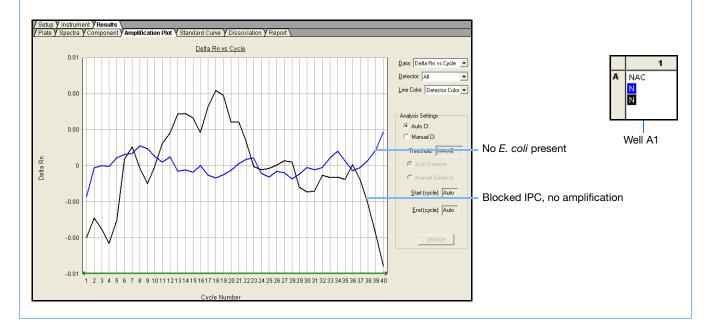


The following Example Experiment boxes show the amplification plots of each type of result (NAC, NTC, plus (+), minus (-), and undetermined (?)).

**Note:** The amplification plots are obtained from the AQ plate document used to amplify the samples in the 96-well plate.

### **Example Experiment**

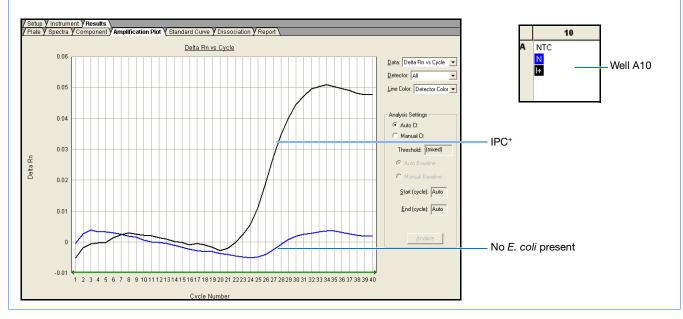
The image below displays the amplification plot for well A1, the No Amplification Control (NAC), which contains blocked IPC and no target template. This plot demonstrates that there is no amplification for IPC or the *E. coli* target sequence.





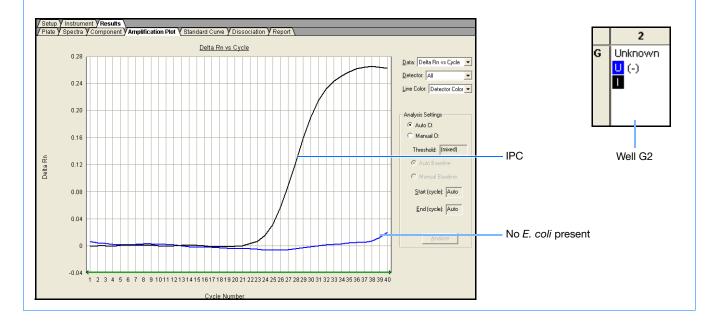
### **Example Experiment**

The image below shows the amplification plot for well A10, the No Template Control (NTC), which contains IPC but no target template. This plot demonstrates a positive amplification curve for IPC and no amplification for the E. coli target sequence.



### **Example Experiment**

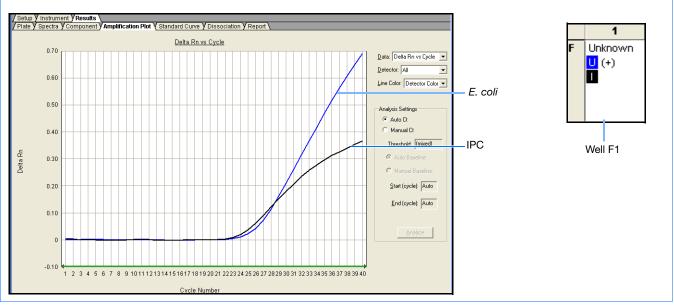
The image below shows the amplification plot for unknown sample well G2, which displays a negative (-) result. The sample is negative because the amplification for the E. coli target sequence is less than the target threshold, while the IPC sample signal is above the calculated IPC threshold defined by the SDS software.





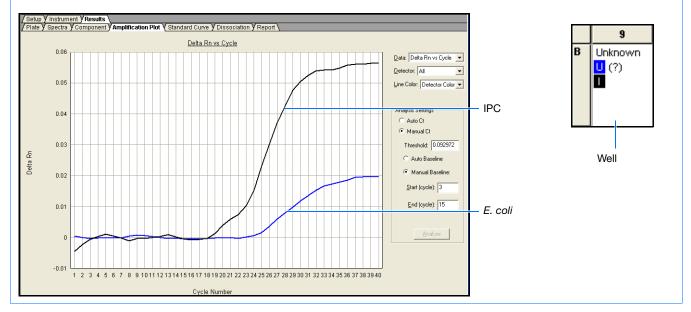
### **Example Experiment**

The image below shows the amplification plot for unknown sample well F1, which displays a positive (+) result. The sample is positive for the *E. coli* target sequence because it has amplification above the target threshold.



### **Example Experiment**

The image below shows the amplification plot for unknown sample well B9, which displays a questionable (?) result. The unknown sample is labeled with a question mark (?) because the unknown sample signal is below the target threshold and the IPC sample signal is below the calculated IPC threshold defined by the SDS software.



**Note:** For more information on analyzing the amplification data and your plus/minus results, see Appendix B, "Viewing Amplification Data," on page 49.

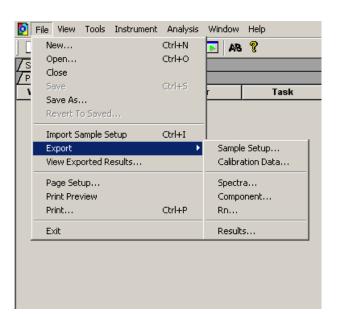
# **Exporting Plus/Minus Plate Data**

You can export numeric data from plus/minus plates into text files, which can then be imported into spreadsheet applications such as Microsoft Excel<sup>®</sup>. You can export graphs as a Microsoft<sup>®</sup> PowerPoint<sup>®</sup> presentation or as JPEG files.

**Note:** You must have PowerPoint installed for the export graphs to PowerPoint feature to work.

- 1. Select **File** > **Export**, then select the data type to export:
  - Sample Setup (\*.txt)
  - Calibration Data (\*.csv)
  - Spectra (\*.csv)
  - Component (\*.csv)
  - Rn (\*.csv)
  - Results (\*.csv)

Refer to the Online Help for information about the export file types.





**2.** Type a file name for the export file.

**Note:** The name of the dialog box depends on the type of data you selected to export.

3. Click Save.

To export data for selected wells and/or report columns to a spreadsheet application:

- 1. Select File > Export > Results.
- **2.** Enter a file name for the export file.
- **3.** Click **Save**. The Export Settings dialog box opens.
- **4.** (Optional) Select export settings:
  - Export only selected wells
  - Apply Report Settings for Data Columns to export the columns selected in the "Report Settings" dialog box (see "Report Tab" on page 55).
- 5. Click OK.

### To export graphs to PowerPoint:

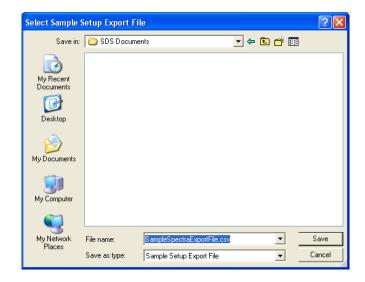
Select Tools > Graph Export > All to
 PowerPoint (or right-click any graph or plate, then select Export All To PowerPoint).

The All to PowerPoint option exports screenshots from all tabs (except the Results > Report tab) of the active file.

Note: To export only the current view, select **Tools** > **Graph Export** > **To PowerPoint** in any view (or right-click any graph or plate, then select **Export To PowerPoint**).

**2.** When prompted, click **OK** to export to PowerPoint. PowerPoint opens and displays your presentation.

**Note:** Title and document information slides are automatically added to your presentation.









- **3.** (Optional) In PowerPoint, modify your presentation.
- **4.** In PowerPoint, click (Save) to save your presentation.

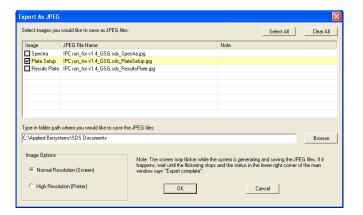
To export plate views or graphs as JPEG files:

1. Select Tools > Graph Export > As JPEG (alternately, right-click any graph or plate, then select Export as JPEG).

The Export as JPEG dialog box opens.

**Note:** In the Export as JPEG dialog box, you can change default file names, select image resolution, and select which plate views or graphs to export, and where the file(s) are saved. Refer to Online Help for more information about this dialog box.

2. Click OK.



# Chapter 6 Performing the Plus/Minus Post-Read Run Exporting Plus/Minus Plate Data

# **Creating Detectors**

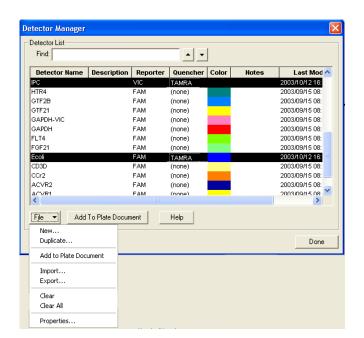
Before you can use a plate document to run a plate, create and apply detectors for all samples on the plate. A detector is a virtual representation of a gene- or allele-specific nucleic acid probe reagent used for analyses performed on instruments.

### To create a detector:

1. Select Tools > Detector Manager.

**Note:** A plate document (any type) must be open before you can access the Tools menu.

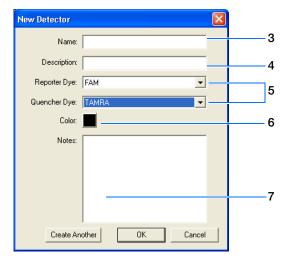
2. Select File > New.



**3.** Enter a name for the detector.

**IMPORTANT!** The name of the detector must be unique and should reflect the target locus of the assay (such as IPC or *E. coli*). Do not use the same name for multiple detectors.

**4.** Optionally, click the **Description** field, then enter a brief description of the detector.



**5.** In the Reporter Dye and Quencher Dye drop-down lists, select the appropriate dyes for the detector.

**Note:** The dyes that appear in the Reporter and Quencher Dye lists are those that have been previously entered using the Dye Manager. If the dye that you want to use does not appear in a list, use the Dye Manager to add the dye and then return to step 5 in this procedure. Refer to the Online Help for more information.

- **6.** Click the **Color** box, select a color to represent the detector using the Color dialog box, then click **OK**.
- **7.** Optionally, click the **Notes** field, then enter any additional comments for the detector.
- **8.** Click **OK** to save the detector and return to the Detector Manager.
- **9.** Repeat steps 2 through 8 for the remaining detectors.
- **10.** In the Detector Manager, click **Done** when you finish adding detectors.

### **Example Experiment**

In the example plus/minus experiment, a detector was created for the *E. coli* target, and another was created for the IPC. The *E. coli* detector was assigned a blue color and IPC a black color.

**Note:** When creating detectors, you use the reporter dye and quencher information (and optionally, the gene name or symbol for the sample name). You can view the contents in a spreadsheet program, such as Microsoft® Excel.

Notes		

# Viewing Amplification Data

# **Configuring Analysis Settings**

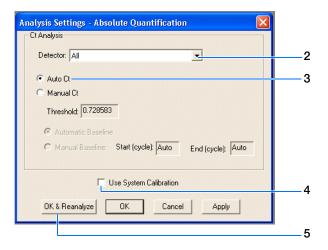
Before you analyze, specify parameters to enable auto-baseline and auto-threshold calculations.

Unless you have already determined the optimal baseline and threshold settings for your experiment, analyze data twice: first using the automatic baseline and threshold feature of the SDS software (Auto  $C_T$ ), and again after determining the optimal baseline and threshold for your data.

### To specify analysis settings:

- 1. Click or select Analysis > Analysis Settings.
- **2.** In the Detector drop-down list, select **All**.
- **3.** Select **Auto Ct** to set the SDS software to automatically generate baseline and threshold values for each well in the study.

**Note:** After analysis, you must verify that the baseline and threshold were called correctly for each well. Alternatively, you can select Manual Ct and specify the threshold and baseline manually. For more information about manually adjusting  $C_T$ , refer to the Online Help.



### Appendix B

Configuring Analysis Settings

**4.** (Optional) Select **Use System Calibration** to use the calibration files that are stored on the computer you are currently using.

**Note:** If you do not select **Use System Calibration**, the calibration information stored in your plate document is used. This information comes from the computer used for data collection when the plate was run.

For more information about system calibration files, refer to the Online Help.

5. Click OK & Reanalyze.

Notes		

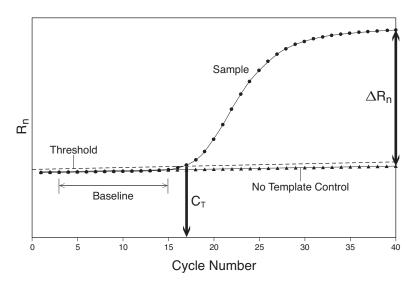
# **Analyzing the Plus/Minus Amplification Data**

Terms Used in Quantitation Analysis

The following are terms commonly used in quantitation analysis.

Term	Definition
Baseline	A line fit to fluorescence intensity values during the initial cycles of PCR, in which there is little change in fluorescence signal.
Threshold cycle (C <sub>T)</sub>	The fractional cycle number at which the fluorescence intensity exceeds the threshold intensity.
Passive reference	A dye that provides an internal fluorescence reference to which the reporter dye signal can be normalized during data analysis.  Normalization is necessary to correct for fluorescence fluctuations caused by changes in concentration or volume.
Reporter dye	The dye attached to the 5' end of a TaqMan® probe. The dye provides a signal that indicates specific amplification.
Normalized reporter (R <sub>n</sub> )	The ratio of the fluorescence intensity of the reporter dye signal to the fluorescence intensity of the passive reference dye.
Delta $R_n (\Delta R_n)$	The magnitude of the signal generated by a set of PCR conditions $(\Delta R_n = R_n - \text{baseline}).$

The figure below is a representative DNA amplification plot and includes some of the terms defined above.



Starting the Analysis

To analyze the amplification data [(AQ) plate], click or select **Analysis > Analyze**. The software generates several types of result views, as described in the following section.

# Viewing the Amplification Data

# About the Results Tab

In the Results tab, you can view the results of the amplification run, change the parameters, and reanalyze the data.

The Results tab has seven secondary tabs. Details about each tab are provided in the Online Help.

- To move between views, click a tab.
- To select all 96 wells on a plate, click the upper-left corner of the plate.
- To adjust graph settings, double-click the y- or x-axis of a plot to display the Graph Settings dialog. The adjustable settings depend on which plot you are viewing.

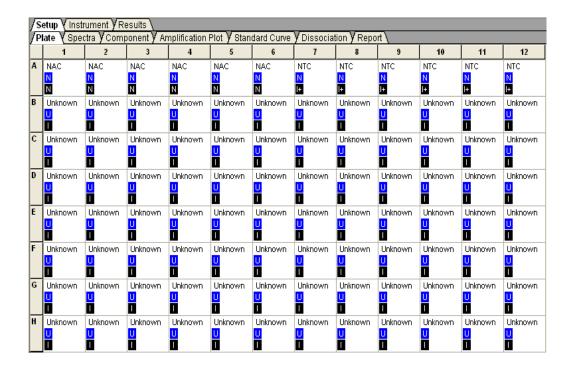


### Plate Tab

Displays the results data of each well, including:

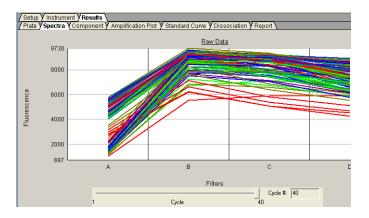
- The sample name and detector task and color for each well.
- A calculated value indicating quantity,  $\Delta Rn$ , or  $C_T$ . *Quantity* is the default. Select **Analysis** > **Display** to select the value to display.

**Note:** For detectors without standards, the Plate Tab displays "Undet." (meaning undetermined).



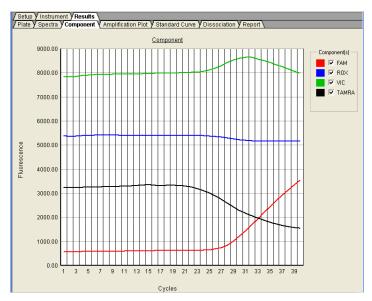
### **Spectra Tab** Displays the fluorescence spectra of selected wells.

- The Cycles slider allows you to see the spectra for each cycle by dragging it with the pointer.
- The Cycle # text box shows the current position of the slider.



### Component Tab

Displays the complete spectral contribution of each dye in a selected wells over the duration of the PCR run. Only the first selected well is shown at one time.

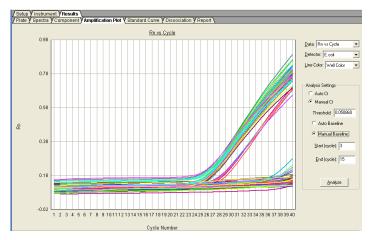


Double-clicking either axis displays the Graph Settings dialog box.

### Amplification Plot Tab

Allows you to view both real-time and post-run amplification of specific samples. The Amplification Plot tab displays all samples in the selected wells.

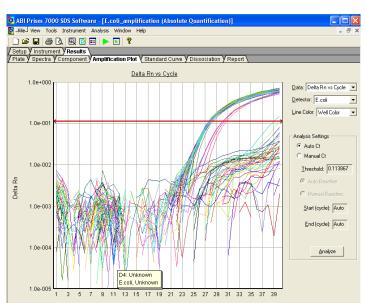
### Rn vs. Cycle (Linear)



This plot displays normalized reporter  $(R_n)$  dye fluorescence as a function of cycle. You can use this plot to identify and examine irregular amplification.

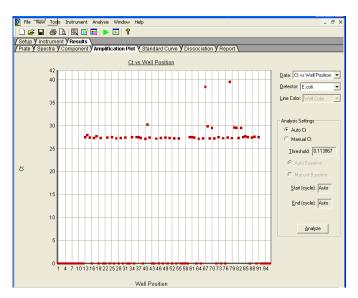
For more information about R<sub>n</sub>, refer to the *Real-Time PCR Systems Chemistry Guide* (PN 4348358).

### △Rn vs.Cycle (Log)



This plot displays the  $R_n$  dye fluorescence as a function of cycle. You can use this plot to identify and examine irregular amplification and to manually set the threshold and baseline parameters for the run.

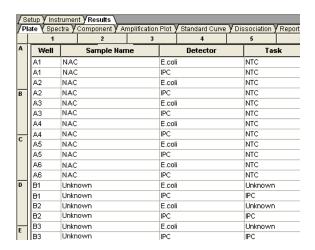
### Ct vs. Well Position Plot



This plot displays threshold cycle  $(C_T)$  as a function of well position. You can use this plot to locate outliers from detector data sets.

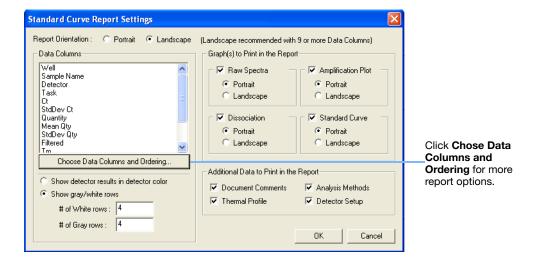
### Report Tab

Displays data for selected wells in a table format. The data columns associated with the report are determined by the assay type. For AQ plate documents, the following data columns are available: Well, Sample Name, Detector, Task, Ct, Std Dev Ct, Quantity, Mean Qty, StdDev Qty, Filtered, Tm, and three User-Defined columns. Refer to Online Help for information on configuring the User-Defined columns.



**Note:** To select the column used to sort the data, click the column heading to sort in ascending (first and alternating clicks) or descending alphanumeric order.

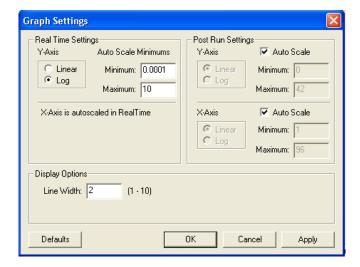
In the Report Settings dialog box, you can format the display of the report and how the report is printed. You have the option (see "Exporting Plus/Minus Plate Data" on page 43) to apply these report settings when you export data. Refer to the Online Help for more information.



### Adjusting Graph Settings

Right-click the Spectra, Component, Amplification Plot, or Standard Curve graph, then select **Graph Settings** to display the Graph Settings dialog box.

The adjustable settings depend on which graph you are viewing. Refer to the Online Help for more information.



# **Example Plus/Minus Experiment**

### Overview

To illustrate how to design, perform, and analyze plus/minus experiments, this section guides you through an example experiment. The example experiment, presented here and in Example Experiment boxes throughout this guide, represents a typical plus/minus experiment setup. You can use the example experiment as a quick-start procedure to familiarize yourself with the plus/minus workflow. See Chapters 1 through 6 for detailed plus/minus workflow procedures.

### Description

The objective of the example plus/minus experiment is to determine if an E. coli target sequence is present or not present in each batch of hamburger meat. The experiment uses duplex PCR where a set of primers and a VIC® dye-labeled probe for the IPC plus a set of primers and a FAM<sup>TM</sup> dye-labeled probe for the target E. coli sequence are run together in each reaction. The set of primers/probe for detecting E. coli was custom designed by Applied Biosystems Primer Express® software.

Reactions were set up for PCR using the TaqMan® Universal PCR Master Mix and appropriate primers and probes.

The example plus/minus experiment data and results were generated using a 7300 system by performing:

- A pre-read run on a plus/minus plate to determine the baseline fluorescence associated with primers and probes before amplification.
- An amplification run using an AQ plate document to generate real-time PCR data, which can be used to analyze and troubleshoot the PCR data for the plus/minus assay, if needed.
- A post-read run using the original plus/minus plate document, which automatically subtracts the pre-read baseline fluorescence determined during the pre-read run, then assigns positive or negative calls using the amplified data.

Notes\_\_\_\_\_

# Example Plus/Minus Experiment Procedure

### Design the experiment and prepare DNA:

- **1.** Design the experiment as explained in Chapter 2.
  - a. Order the TaqMan® Exogenous Internal Positive Control Reagents kit and the TaqMan® Universal PCR Master Mix.
  - b. Design the primers and FAM<sup>™</sup> dye-labeled probe set for *E. coli* detection with Applied Biosystems Primer Express software.
- 2. Extract DNA from samples (see "Preparing DNA" on page 12) using the PrepMan® Ultra Sample Preparation Reagent with Protocol (PN 4322547) to obtain a final concentration of 10 ng/μL of DNA for each sample.
- **3.** Prepare sufficient reaction mix (see "Preparing the PCR Reaction Mix" on page 14) by using the volumes as listed in the table on the right

CAUTION CHEMICAL HAZARD.

TaqMan® Universal PCR Master Mix (2×) No AmpErase® UNG may cause eye and skin irritation. Exposure may cause discomfort if swallowed or inhaled. Read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Item	Volume for one Reaction (Standard plate)	Volume for one Reaction (Fast plate)	Final Concentration
TaqMan <sup>®</sup> Universal PCR Master Mix (2X)	25.0	10.0	1×
10X Exo IPC Mix (IPC kit)	5.0	2.0	50 to 900 nM
50X Exo IPC DNA (IPC kit)	1.0	0.4	50 to 900 nM
Target primers, probe, and deionized water	14.0	5.6	50 to 250 nM
Total	45.0	18.0	_

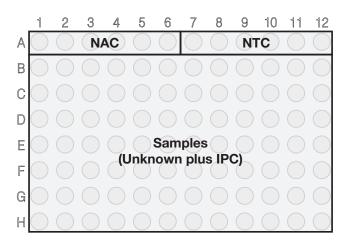
Notes		

### **4.** Prepare the reaction plate:

- a. Pipette 45  $\mu$ L (18  $\mu$ L for the Fast system) of the reaction mixture into each well of a 96-well reaction plate.
- b. Pipette 5 μL (2 μL for the Fast system) of IPC block, TE or water, or unknown sample into the designated wells of a 96-well plate such as the example indicated in the table to the right. (see "Setting Up the Reaction Plate" on page 13).

**Note:** The final reaction volume in each well is  $50 \,\mu\text{L}$  (20  $\mu\text{L}$  for the Fast system).

**c.** Keep the reaction plate on ice until you are ready to load it into the 7300/7500/7500 Fast system.



Wells	If preparing	Add (Standard plate)	Add (Fast plate)
A1 to A6	NAC <sup>‡</sup>	5 μL of 10X Exo IPC Block	2 μL of 10X Exo IPC Block
A7 to A12	NTC <sup>§</sup>	5 μL of 1X TE or H <sub>2</sub> O	2 μL of 1X TE or H <sub>2</sub> O
B1 to H12	U#	5 μL of sample being tested for <i>E. coli</i>	2 μL of sample being tested for E. coli

- ‡ No Amplification Control Well contains no target template and no IPC.
- § No Template Control Well contains no target template, only IPC.
- # Unknown Well contains both target template and IPC.

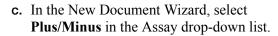
00	-	. •	, c

### Perform the pre-read run:

**1.** Create a plus/minus plate document:

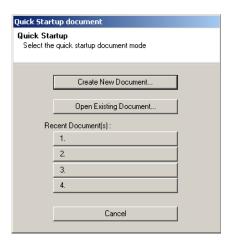
Follow the instructions as described in Chapter 4. Briefly,

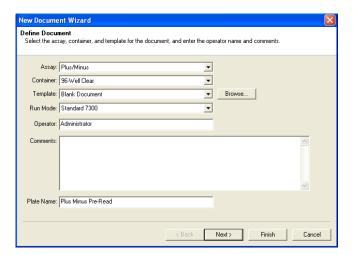
- a. Select Start > All Programs > Applied Biosystems > 7300/7500/7500 Fast System > 7300/7500/7500 Fast System Software () to start the 7300/7500/7500 Fast SDS software.
- **b.** In the Quick Startup document dialog box, select **Create New Document**.



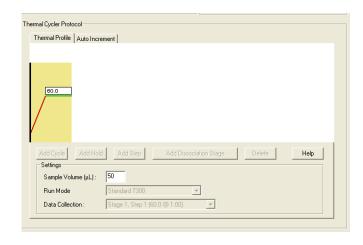
- d. In the Plate Name field, type Plus Minus Pre-Read, then click Next.
- e. Add detectors to the plate document (see Appendix A, Creating Detectors on page 47) then click **Next**.
- **f.** Specify detectors and tasks for each well, then click **Finish**.
- **g.** Double-click each well to type the sample name, then save the document.
- 2. Enter the sample names and specify tasks in the Well Inspector (View > Well Inspector).

**IMPORTANT!** If your experiment does not use all the wells on a plate, do not omit the wells from use at this point. You can omit unused wells after the run is completed. For more information about omitting wells, refer to the Online Help.





- **3.** Perform the plus/minus pre-read run.
  - **a.** Select the **Instrument** tab. By default, the standard PCR conditions are displayed.
  - b. Select **File > Save**, type a name for the plus/minus plate document, then click **Save**.
  - **c.** Load the reaction plate into the instrument.
  - d. Click Pre-Read.



## Amplify the DNA:

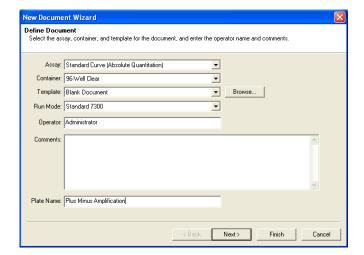
**1.** Create an AQ plate document for amplification:

Follow the instructions as described in Chapter 5. Briefly, in the New Document Wizard,

- a. Select Standard Curve (Absolute Quantitation) in the Assay drop-down list.
- b. In the Plate Name field, type Plus Minus Amplification, then click Next.

**Note:** A standard curve is not needed for a non-quantitation amplification run.

- **c.** Add detectors to the plate document (see Appendix A, "Creating Detectors," on page 47), then click **Next**.
- **d.** Specify the detectors and tasks for each well, then click **Finish**.
- **e.** Type the sample names, then save the document.

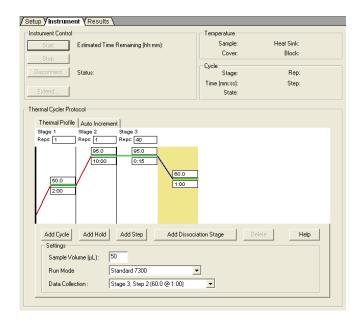


Notes

- **2.** Perform the plus/minus amplification run.
  - a. Select the **Instrument** tab.
  - **b.** Select **File > Save**, type a name for the AQ plate document, then click **Save**.
  - **c.** Load the reaction plate into the instrument, then click **Start**.

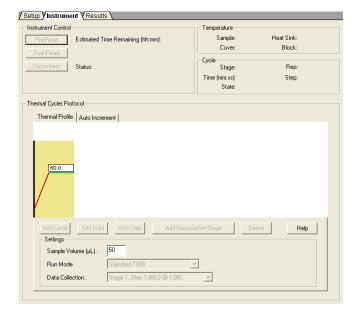
By default, the standard PCR conditions for the PCR step are displayed.

After the run, a message indicates if the run is successful, or if errors were encountered.



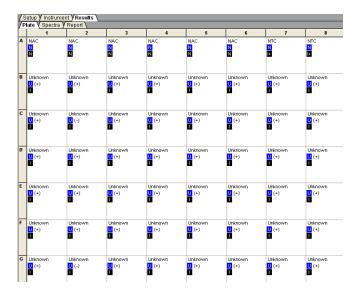
### Perform the plus/minus post-read run:

- 1. Open the plus/minus pre-read plate document, then use it to perform the post-read run. (See "Performing the Post-Read Run" on page 36.)
  - a. Select the **Instrument** tab.
  - b. Select File > Save As, type a name for the plus/minus post-read plate document, then click Save.
  - **c.** Load the reaction plate into the instrument.
  - d. Click Post-Read.
  - e. Click or select Analysis > Analyze. Click the **Results** tab to view results for each well.



Notes

**2.** If you need to troubleshoot the plus/minus results, see "Viewing the Amplification Data" on page 52.



Notes



## References

Kwok, S. and Higuchi, R. 1989. Avoiding false positives with PCR. *Nature* 339:237–238.

Mullis, K.B. and Faloona, F.A. 1987. Specific synthesis of DNA *in vitro* via a polymerase-catalyzed chain reaction. *Methods Enzymol.* 155:335–350.

Saiki, R.K., Scharf, S., Faloona, F., *et al.* 1985. Enzymatic amplification of  $\beta$ -globin genomic sequences and restriction site analysis for diagnosis of sickle cell anemia. *Science* 230:1350–1354.

References

# Index

Numerics	hazards xiv
7300/7500 system	chemical waste disposal, guidelines xiv
description 2	Color box 48
plus/minus assay 2	Component tab 53
7300/7500/7500 Fast system	conventions
background and pure-dye runs 18	bold text vii
96-well reaction plate 13, 59 loading 25	for describing menu commands vii IMPORTANTS! vii in this guide vii italic text vii
	Notes vii
absolute quantitation 61	user attention words vii
amplification data	Ct vs. Well Position view 55
generating 27	Ct. See threshold cycle
viewing 52 amplification plates detectors, creating 47	customer feedback, on Applied Biosystems documents viii
starting a run 34	D
amplification plot 40, 41	_
Amplification Plot tab 54	DANGER, description xi
amplification run	delta Rn 51
PCR conditions 33	Delta Rn vs. Cycle view 54
performing 33	Detector Manager window 47
Analysis button 34	detector tasks 19
analysis settings 49	detectors 18, 61
appearance of graphs 52	amplification run 30
Applied Biosystems	creating 47 definition 47
customer feedback on documentation viii	in document window 21
Information Development department viii	removing 30
assumptions for using this guide vii auto Ct 49	selecting 21
	display options 52
autoscaling options 52	DNA
В	amplifying 51, 61 chemistry systems 12
baseline 49, 51	final concentration 12 isolation 12, 58
biological hazard guidelines xiii	quality 12
bold text, when to use vii	documentation, related viii
	documents, plate
C	importing 20
calibration, background and pure dye 50	plus/minus plate document 18 templates 20
CAUTION, description vii, xi chemical safety guidelines xii	dyes 51

chemical waste

E	N
exporting plate data 43	New Detector window 47
	New Document Wizard window 20, 60
F	no amplification control 21, 30
FAM-labeled probe 9, 58	See also NAC 3, 16, 18 plot 40
fluorescence	no template control 22, 31
spectra 18, 53	See also NTC 3, 16, 18
fluorogenic	normalized reporter 51
5' nuclease chemistry 8	Notes field 48
G	nucleotide sequence 3
graph settings 52	
Graph Settings dialog box 56	0
guidelines	OK & Reanalyze button 50
chemical safety xii	omitting unused wells 23
chemical waste disposal xiv	online help 60
	options, display 52
H	options, graph 52
hazard icons xi	
hazards	P
chemical waste xiv	passive reference 31, 51
	(ROX dye) 23, 32
1	PCR
IMPORTANT, description xi	cycles 8
Information Development department, contacting viii	duplex 57
Instrument tab 24, 33, 36, 61, 62	reaction mix 14
IPC	real-time 28, 34 starting an amplification plate run 34
amplification plot 41	plate document
blocked 21, 30	amplification 28
blocking reagent 59	creating 18
definition 2	creating for AQ 29
TaqMan Exogenous Internal Positive Control Reagents Kit 2	parameters 18
VIC probe 2	sample names 22 Plate tab 38, 39, 52
VIC reporter dye 21	
italic text, when to use vii	plates, standard vs. fast 15
	plot appearance 52
L	plus/minus assay about 4
line width 52	amplification data 51
	amplification run 61
M	calling unknowns using IPC 39
	calls 18
manual Ct 49	chemistry kits 9 controls and unknowns 18, 59
menu commands, conventions for describing vii	end-point assay 28
MSDSs	final reaction volume 16
referring to xiii	IPC 2
MSDSs, obtaining ix	negative result 41
	plate document 18, 19, 60 post-read run 36, 62
	probe and primers 9

reaction mix 58 reaction plate 59	settings, graph 52
result calls 39, 62	Setup tab 23 software 18
sample experiment 57	
tasks 19 terms used 3	Spectra tab 53
using absolute quantitation for real-time data 29	starting an amplification plate run 34
viewing results 39	<b>-</b>
workflow 4	Т
plus/minus plate documents	TAMRA dye-labeled quencher 9
exporting plus/minus data 43 importing sample information 20 templates 20	TaqMan Exogenous Internal Positive Control Reagents Kit 9, 58
Post-Read button 37, 62	TaqMan probe-based chemistry 8
post-read run	TaqMan Universal PCR Master Mix 9, 13, 57, 58
analysis 62	protocol 9
performing 36	target template 3, 21
PrepMan Ultra Sample Preparation Reagent Kit 12, 58	task 18
Pre-Read button 25, 61	template documents 20
pre-read run 25, 60	text conventions vii
creating plate document 18, 20	threshold 49
performing 24	threshold cycle
Primer Express software 9, 58	auto Ct 49
probe and primers, designing 9	definition 51 Tools menu 47
Q	training, information on ix
quantitation analysis, terms 51	troubleshooting 28, 63
Quencher Dye drop-down list 48	U
R	unknown sample 3, 22
	unknowns 16
reaction plate 59 setting up 13	Use checkbox 30
setting up 13 real-time PCR 28	Use System Calibration checkbox 50
Real-Time PCR Systems Chemistry Guide 2	user attention words, described vii
reference, passive 31, 51	
Report tab 55	V
reporter dye 51	VIC-labeled probe 9
Reporter Dye drop-down list 48	viewing results 39, 52
required materials 4	viewing results 37, 32
•	W
Results tab 34, 52, 62	
Rn vs. Cycle view 54	WARNING, description vii, xi
Rn. See normalized reporter	Well Inspector window 22, 60
ROX dye 31	width, lines 52
S	X
safety	x-axis 52
biological hazards xiii chemical waste xiv	Υ
sample experiment 5, 9, 12, 13, 16, 23, 38, 40, 41, 42,	y-axis 52
48, 57, 58	y-anis 32

Index

### **Worldwide Sales and Support**

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

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

### Headquarters

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

06/2010

