

***ABI PRISM<sup>®</sup>  
BigDye<sup>™</sup> Primer  
v3.0 Ready Reaction  
Cycle Sequencing  
Kit***

*For –21 M13 and M13 REV Primers*

---

---

**Protocol**

---

---

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

**AB** Applied  
Biosystems

© Copyright 2001, 2010 Applied Biosystems

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

**Notice to Purchaser: Limited License**

The purchase of the ABI PRISM® BigDye™ Primer v3.0 Ready Reaction Cycle Sequencing Kit includes a limited, nontransferable, non-exclusive license (without the right to resell, repackage, or sublicense) under U.S. patents and corresponding foreign patents and patent applications to use this product solely with an Applied Biosystems commercial automated DNA sequencing machine or other authorized automated DNA sequencing machines that have been authorized under this patent by Applied Biosystems. No license is hereby granted for the use of this kit, or the reagents therein, in any other automated sequencing machine. Such license is granted solely for research and other uses that are not unlawful. No other license is granted expressly, impliedly, or by estoppel. For information concerning the availability of additional licenses to practice the patented methodologies, contact: Director of Licensing at Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404. Patents are pending in countries outside the United States.

**Notice to Purchaser: Limited License**

The purchase price of this product includes a limited, nontransferable license under U.S. patents or its foreign counterparts, to use only this amount of the product for DNA Sequencing and related processes described in said patent solely for the research and development activities of the purchaser. No license under these patents to use the PCR process is conveyed expressly or by implication to the purchaser by the purchase of this product. A license to use the PCR Process for certain research and development activities accompanies the purchase of certain reagents from licensed suppliers such as Applied Biosystems when used in conjunction with an Authorized Thermal Cycler, or is available from Applied Biosystems. Further information on purchasing licenses to practice PCR Process may be obtained by contacting the Director of Licensing at Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404.

**Notice to Purchaser About Limited License**

This kit (reagent) is sold pursuant to a limited sublicense from Amersham International plc under one or more U.S. Patent Nos. 5,498,523, 5,614,365 and corresponding foreign patents and patent applications. The purchase of this kit (reagent) includes a limited non-exclusive sublicense (without the right to resell, repackage, or further sublicense) under such patent rights to use this reagent for DNA sequencing or fragment length analysis solely with an Applied Biosystems commercial automated sequencing machine or other authorized DNA sequencing machines that have been authorized for such use by Applied Biosystems, or for manual DNA sequencing. No license is hereby granted for the use of this kit, or the reagents therein, in any other automated sequencing machine. Such sublicense is granted solely for research and other uses that are not unlawful. No other license is granted expressly, impliedly, or by estoppel. For information concerning the availability of additional license to practice the patented methodologies, contact: Amersham Life Science, Inc., Vice President, Regulatory Affairs, P.O. Box 22400, Cleveland, Ohio 44122. Patents are pending in countries outside the United States.

ABI PRISM and its design, Applied Biosystems, BioCAD, BioLIMS, BioMerge, Cytofluor, GeneScan, MicroAmp, POROS, Primer Island, and Procise are registered trademarks and AB (Design), Applera, BigDye, Biospectrometry, Catalyst, Expedite, FMAT, Hi-Di, Lane Guide, Mariner, POP-4, POP-5, POP-6, SQL\*GT, and Voyager are trademarks of Applied Biosystems or its subsidiaries in the U.S. and certain other countries.

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

Centricon is a registered trademark of W. R. Grace and Co.

Macintosh is a registered trademark of Apple Computer, Inc.

pGEM is a registered trademark of Promega Corporation.

Long Ranger is a trademark of FMC Corporation.

Windows NT is a registered trademarks of the Microsoft Corporation.

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

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

# Contents

## **1 Introduction**

Chapter Summary .....	1-1
In This Chapter .....	1-1
About the Kit .....	1-2
Reagent Requirements .....	1-2
BigDye Primer v3.0 Ready Reaction Kit .....	1-2
Cycle Sequencing with AmpliTaq DNA Polymerase, FS .....	1-3
BigDye Primer v3.0 Appearance on the 377 or 373 Instrument Gel Image .....	1-3
Instruments .....	1-4
Instrument Platforms .....	1-4
Thermal Cyclers .....	1-4
Required Software .....	1-5
Dye/Filter Sets and Matrix Standards for the 310, 377, and 373 Instruments .....	1-5
Dye Sets and Spectral Standards for the 3700 and 3100 Instruments	1-5
Instructions .....	1-6
Dye Set/Primer (Mobility) Files .....	1-6
Reagents and Storage .....	1-8
Available Kits .....	1-8
Storage and Use of the Kit .....	1-8
Materials Supplied by the User .....	1-9
Overview .....	1-9
Materials for Cycle Sequencing .....	1-10
Materials for Purifying Extension Products .....	1-11
Safety .....	1-11
Documentation User Attention Words .....	1-11

Chemical Hazard Warning . . . . .	1-11
Chemical Waste Hazard Warning . . . . .	1-12
Site Preparation and Safety Guide . . . . .	1-12
About MSDSs . . . . .	1-12
Ordering MSDSs . . . . .	1-13

## ***2 Preparing the Templates***

Chapter Summary . . . . .	2-1
In This Chapter. . . . .	2-1
Control DNA Templates . . . . .	2-2
Using Control DNA . . . . .	2-2
Control DNA Sequence . . . . .	2-2
An Additional Control Sold Separately . . . . .	2-2
Template Preparation Methods . . . . .	2-3
Single- and Double-Stranded Templates . . . . .	2-3
BAC DNA Templates . . . . .	2-3
PCR Templates . . . . .	2-4
Direct Link PCR Sequencing . . . . .	2-4
Purifying PCR Fragments . . . . .	2-5
DNA Quantity . . . . .	2-6
Quantitating DNA . . . . .	2-6
Template Quantity . . . . .	2-6

## ***3 Performing Cycle Sequencing***

Chapter Summary . . . . .	3-1
In This Chapter. . . . .	3-1
Introduction . . . . .	3-2
Overview . . . . .	3-2
Optimizing Cycle Parameters . . . . .	3-2
Cycle Sequencing Single- and Double-Stranded DNA . . . . .	3-3
Overview . . . . .	3-3
Preparing the Reactions . . . . .	3-3
Cycle Sequencing on the System 9700, 9600, or 2400 . . . . .	3-6

Cycle Sequencing on the TC1 or 480 . . . . .	3-7
Cycle Sequencing BAC DNA . . . . .	3-8
Preparing Sequencing Reactions . . . . .	3-8
Performing Cycle Sequencing . . . . .	3-8

#### ***4 Preparing the Samples for Loading***

Chapter Summary . . . . .	4-1
In This Chapter . . . . .	4-1
Recommendations for Precipitation/Loading . . . . .	4-2
For the 3700, 3100, 310, and 373 Instruments . . . . .	4-2
For the 377 Instrument . . . . .	4-2
Ethanol Precipitation for Microcentrifuge Tubes . . . . .	4-3
Procedure . . . . .	4-3
Procedural Options . . . . .	4-4
Ethanol Precipitation for 96-Well Reaction Plates . . . . .	4-5
Standard Method . . . . .	4-5
Precipitating in 96-Well Reaction Plates . . . . .	4-5

#### ***5 Sample Electrophoresis***

Chapter Summary . . . . .	5-1
In This Chapter . . . . .	5-1
Before You Begin . . . . .	5-2
Important Reminders . . . . .	5-2
Electrophoresis on the ABI PRISM 3700 DNA Analyzer . . . . .	5-3
Requirements . . . . .	5-3
Performing Sample Electrophoresis . . . . .	5-3
Electrophoresis on the ABI PRISM 3100 Genetic Analyzer . . . . .	5-4
Requirements . . . . .	5-4
Performing Sample Electrophoresis . . . . .	5-4
Electrophoresis on the ABI PRISM 310 Genetic Analyzer . . . . .	5-5
Requirements . . . . .	5-5
Setting the Data Analysis Range . . . . .	5-6
Resuspending the Samples for Sequencing with POP-6 Polymer . . . . .	5-6

Resuspending the Samples for Sequencing with POP-4 Polymer . . .	5-7
Electrophoresis on the ABI PRISM 377 DNA Sequencers . . . . .	5-8
Requirements . . . . .	5-8
Using the Lane Guide Kit . . . . .	5-9
Using Long-Read Gel and Buffer Formulations . . . . .	5-9
Resuspending and Loading the Samples . . . . .	5-10
Electrophoresis on the 373 Instrument with BigDye Filter Wheel. . . . .	5-12
Requirements . . . . .	5-12
Resuspending and Loading the Samples . . . . .	5-13

## ***A Control DNA Sequences***

Appendix Summary . . . . .	A-1
In This Appendix . . . . .	A-1
-21 M13 Forward Primer . . . . .	A-2
Sequence . . . . .	A-2
M13 Reverse Primer . . . . .	A-3
Sequence . . . . .	A-3

## ***B Technical Support***

Technical Support . . . . .	B-1
Contacting Technical Support . . . . .	B-1
To Contact Technical Support by E-Mail. . . . .	B-1
Hours for Telephone Technical Support . . . . .	B-2
To Contact Technical Support by Telephone or Fax . . . . .	B-2
To Reach Technical Support Through the Internet . . . . .	B-6
To Obtain Documents on Demand. . . . .	B-7

# *Introduction*

# 1

## Chapter Summary

---

**In This Chapter** The following topics are covered in this chapter:

<b>Topic</b>	<b>See Page</b>
About the Kit	1-2
Instruments	1-4
Required Software	1-5
Reagents and Storage	1-8
Materials Supplied by the User	1-9
Safety	1-11

---

## About the Kit

---

**Reagent Requirements**

- The ABI PRISM® BigDye™ Primer v3.0 Ready Reaction Cycle Sequencing Kit with AmpliTaq® DNA Polymerase, FS, requires **new** instrument (matrix) files for the ABI PRISM® 310 Genetic Analyzer, ABI PRISM® 377 DNA Sequencers, and ABI PRISM® 373 DNA Sequencers with the ABI PRISM® BigDye™ Filter Wheel installed and **new** spectral calibrations for the ABI PRISM® 3700 DNA Analyzer and the ABI PRISM® 3100 Genetic Analyzer:
  - The 310, 377, and 373 instruments require the ABI PRISM® BigDye™ Matrix Standards v3.0 (P/N 4390421) for instrument (matrix) file generation.
  - The 3700 instrument requires ABI PRISM® 3700 BigDye™ Terminator v3.0 Sequencing Standard (P/N 4390309) for spectral calibration.
  - The 3100 instrument requires ABI PRISM® BigDye™ Terminator v3.0 Sequencing Standard (P/N 4390303) for spectral calibration.
- The dRhodamine Matrix Standards and Matrix Standard Set DS-01 are not compatible with BigDye™ primers v3.0, BigDye™ terminators v3.0, or dGTP BigDye™ terminators v3.0.
- The alcohol precipitation methods are different from those recommended for the original BigDye primers.
- There are new v3.0 mobility files.
- The basecallers are the same.

---

**BigDye Primer v3.0 Ready Reaction Kit**

The ABI PRISM BigDye Primer v3.0 Ready Reaction Cycle Sequencing Kit combines AmpliTaq® DNA Polymerase, FS, the BigDye™ primers v3.0, and all the required components for the sequencing reaction.

The kit contains nucleotide mixes that have been specifically optimized for AmpliTaq DNA Polymerase, FS. The deoxy- and dideoxynucleotide ratios in the nucleotide mixes have been formulated to give a good signal balance above 700 bases. These formulations also contain 7-deaza-dGTP in place of dGTP to minimize band compressions.

In the Ready Reaction format, the dye-labeled primers, deoxynucleoside triphosphates, dideoxynucleoside triphosphates, AmpliTaq DNA Polymerase, FS, and buffer are premixed into A, C, G,



and T Ready Reaction cocktails to eliminate time-consuming reagent preparation. These reagents are suitable for performing fluorescence-based cycle sequencing reactions on single-stranded or double-stranded DNA templates, on polymerase chain reaction (PCR) fragments, and on large templates, *e.g.*, the ends of BAC clones.

---

---

**Cycle Sequencing  
with  
AmpliTaq DNA  
Polymerase, FS**

The kit formulation contains the sequencing enzyme AmpliTaq DNA Polymerase, FS. This enzyme is a variant of *Thermus aquaticus* DNA polymerase that contains a point mutation in the active site. This results in less discrimination against dideoxynucleotides and a much more even peak intensity pattern.

This enzyme also has a second mutation in the amino terminal domain that virtually eliminates the 5'→3' nuclease activity of AmpliTaq DNA Polymerase. The enzyme has been formulated with a thermally stable inorganic pyrophosphatase to eliminate problems associated with pyrophosphorolysis.

Cycle sequencing protocols that rely on the use of AmpliTaq DNA Polymerase, FS offer the following advantages over traditional sequencing methods:

- Less hands-on operation
- No alkaline denaturation step required for double-stranded DNA
- Same protocol for both single- and double-stranded templates
- Less starting template needed
- More reproducible results

---

---

**BigDye Primer  
v3.0 Appearance  
on the 377 or 373  
Instrument Gel  
Image**

The dye/base relationships and colors of the BigDye primers v3.0 as they appear on the gel image are shown below for the 377 and 373 instruments.

Base	Primer	Color of Bands on ABI PRISM 377 or 373 Instrument Gel Image
A	V3 Dye 2	Green
C	V3 Dye 1	Blue
G	V3 Dye 3	Yellow
T	V3 Dye 4	Red

## Instruments

---

**Instrument Platforms** The ABI PRISM BigDye Primer v3.0 Ready Reaction Cycle Sequencing Kit is for use with the following instruments:

- ABI PRISM® 3700 DNA Analyzer
- ABI PRISM® 3100 Genetic Analyzer
- ABI PRISM® 310 Genetic Analyzer
- ABI PRISM® 377 DNA Sequencer (all models<sup>1</sup>)

This kit can also be used with ABI PRISM® 373 DNA Sequencers with the new ABI PRISM® BigDye Filter Wheel installed.<sup>2</sup> Refer to the *ABI PRISM BigDye Filter Wheel User Bulletin* (P/N 4304367) for more information.

General instructions are given for using the kit reagents to generate samples for these instruments. For more detailed instructions, refer to the appropriate instrument user's manual.

**IMPORTANT** This kit is not designed for use with ABI PRISM® 373 DNA Sequencers and ABI PRISM® 373 DNA Sequencers with XL Upgrade that do not have the ABI PRISM BigDye Filter Wheel.

---

**Thermal Cyclers** The protocols provided in this document were optimized using Applied Biosystems thermal cyclers, including:

- GeneAmp® PCR Instrument Systems 9700, 9600, and 2400
- DNA Thermal Cycler 480
- DNA Thermal Cycler (TC1)

If you use a thermal cycler not manufactured by Applied Biosystems, you may need to optimize thermal cycling conditions. Ramping time is very important. If the thermal ramping time is too fast (>1°/sec), poor (noisy) data may result.

---

1. Includes the ABI PRISM 377, ABI PRISM 377-18, ABI PRISM 377 with XL Upgrade, and the ABI PRISM 377 with 96-Lane Upgrade instruments.

2. Includes the ABI PRISM 373 and ABI PRISM 373 with XL Upgrade instruments.

## Required Software

**Dye/Filter Sets and Matrix Standards for the 310, 377, and 373 Instruments** The dye/filter sets and matrix standards required for the 310, 377, and 373 instruments are listed in the table below.

**IMPORTANT** The instrument (matrix) file for the BigDye primers v3.0 cannot be used for the BigDye terminators (original), BigDye terminators v2.0, dRhodamine terminators, or BigDye primers (original).

Instrument	Dye/Filter Set	Standards for Instrument (Matrix) File Generation
310 Genetic Analyzer	Filter Set E	ABI PRISM® BigDye™ Matrix Standards v3.0 (P/N 4390421)
377 DNA Sequencers <sup>a</sup>	Filter Set E	
373 DNA Sequencers with the BigDye Filter Wheel <sup>b</sup>	Filter Set A	

a. Includes the ABI PRISM 377, ABI PRISM 377-18, ABI PRISM 377 with XL Upgrade, and the ABI PRISM 377 with 96-Lane Upgrade instruments.

b. Includes the ABI PRISM 373 and ABI PRISM 373 with XL Upgrade instruments.

**Dye Sets and Spectral Standards for the 3700 and 3100 Instruments** The dye sets and spectral standards required for the 3700 and 3100 instruments are listed in the table below.

**IMPORTANT** Spectral calibrations for the BigDye primers v3.0 are not compatible with the BigDye terminators (original), BigDye terminators v2.0, dRhodamine terminators, or BigDye primers (original).

Instrument	Dye Set	Standards for Spectral Calibration
3700 DNA Analyzer	D	ABI PRISM® 3700 BigDye™ Terminator v3.0 Sequencing Standard (P/N 4390309)
3100 Genetic Analyzer	Z	ABI PRISM® BigDye™ Terminator v3.0 Sequencing Standard (P/N 4390303)

---

---

**Instructions For Generating Matrices**

- For the 377 and 310 instruments, refer to the product insert for instructions on using the ABI PRISM BigDye Matrix Standards v3.0 (P/N 4390421) to generate matrices.
- For the 373 instruments, contact Technical Support for instructions on using the ABI PRISM BigDye Matrix Standards v3.0 (P/N 4390421) to generate matrices.

**For Performing Spectral Calibrations**

- For the 3700 instrument, refer to the product insert for instructions on using the ABI PRISM 3700 BigDye Terminator v3.0 Sequencing Standard (P/N 4390309) to perform spectral calibration.
- For the 3100 instrument, refer to the product insert for instructions on using the ABI PRISM BigDye Terminator v3.0 Sequencing Standard (P/N 4390303) to perform spectral calibration.

---

---

**Dye Set/Primer (Mobility) Files Available in Two Places**

To analyze sequencing data generated with BigDye chemistries v3.0, you need dye set/primer (mobility) files that were created for v3.0 chemistries. The dye set/primer (mobility) files can be obtained from two places:

- The files can be installed from the two CD-ROMs or one floppy disk enclosed in the v3.0 matrix and sequencing standards listed below. See page 1-7 for instructions.
  - ABI PRISM BigDye Matrix Standards v3.0 (P/N 4390421)
  - ABI PRISM BigDye Terminator v3.0 Sequencing Standard (P/N 4390303)
  - ABI PRISM 3700 BigDye Terminator v3.0 Sequencing Standard (P/N 4390309)
- The files can be downloaded from the Internet. See page 1-7 for instructions.

### **Installing Files from the CD-ROMs or Floppy Disk Enclosed in the v3.0 Matrix and Sequencing Standards**

<b>If you wish to analyze your data using a ...</b>	<b>and your data was collected on a...</b>	<b>Refer to the CD-ROM or floppy disk labeled...</b>
computer with the Windows NT® platform	3700, 3100, 310, or 377 instrument	PN 4326478, For Windows NT platform
Macintosh® computer with a CD-ROM drive	310, 377, 373 instrument	PN 4326479, For Macintosh platform
Macintosh computer with a floppy drive	310, 377, 373 instrument	PN 4326480, For Macintosh platform

### **Downloading Files from the Internet**

Dye set/primer (mobility) files can be downloaded from our website:

<http://www.appliedbiosystems.com/techsupp/swpps/SAsw.html>

If you do not have access to the Internet, you can get the files from Applied Biosystems Technical Support, or from your local field applications specialist (call your local sales office for more information).

---

## Reagents and Storage

---

**Available Kits** The following kits are available:

Kit	Primer	Number of Reactions	Part Number
The ABI PRISM® BigDye™ Primer v3.0 Ready Reaction Cycle Sequencing Kit with AmpliTaq® DNA Polymerase, FS	-21 M13	100	4390161
	M13 reverse	100	4390163
	-21 M13	5000	4390157
	M13 reverse	5000	4390159

---

### Storage and Use of the Kit

- Store the kit at –15 to –25 °C.
- Avoid excess (*i.e.*, no more than 5–10) freeze-thaw cycles. Aliquot reagents in smaller amounts if necessary.
- Before each use of the kit, allow the frozen stocks to thaw at room temperature (do not heat).

**IMPORTANT** Mix each stock thoroughly and then centrifuge briefly to collect all the liquid at the bottom of each tube.

- Whenever possible, thawed materials should be kept on ice during use. Do not leave reagents at room temperature for extended periods.
-

## Materials Supplied by the User

---

**Overview** In addition to the reagents supplied in this kit, other items are required.

This section lists general materials needed for:

- Cycle sequencing
- Purifying extension products

**Note** Many of the items listed in this section are available from major laboratory suppliers (MLS) unless otherwise noted. Equivalent sources may be acceptable where noted.

Refer to the individual instrument protocols for the specific items needed for each instrument.

**⚠ WARNING CHEMICAL HAZARD.** Before handling the chemical reagents needed for cycle sequencing, read the safety warnings on the reagent bottles and in the manufacturers' Material Safety Data Sheets (MSDSs). Always wear protective equipment (lab coat, safety glasses, and chemical-resistant gloves) when handling chemicals. Dispose of waste in accordance with all local, state/provincial, and national environmental and health regulations.

---

**Materials for Cycle Sequencing** The table below lists the plates or tubes required for the recommended Applied Biosystems thermal cyclers (page 1-4).

<b>Thermal Cycler</b>	<b>Plate or Tube</b>	<b>Applied Biosystems Part Number</b>
GeneAmp® PCR System 9700	MicroAmp® 96-Well Reaction Plate	N801-0560
	MicroAmp® Reaction Tubes, 0.2-mL	N801-0533
	MicroAmp® Caps, 12 or 8/strip	N801-0534 or N801-0535
	ABI PRISM™ Optical Adhesive Cover Starter Pack or ABI PRISM® Optical Adhesive Covers	4313663 or 4311971
GeneAmp® PCR System 9600	MicroAmp® 96-Well Reaction Plate	N801-0560
	MicroAmp® Reaction Tubes, 0.2-mL	N801-0533
	MicroAmp® Caps, 12 or 8/strip	N801-0534 N801-0535
	ABI PRISM™ Optical Adhesive Cover Starter Pack or ABI PRISM® Optical Adhesive Covers	4313663 or 4311971
GeneAmp® PCR System 2400	MicroAmp® Reaction Tubes, 0.2-mL	N801-0533
	MicroAmp® Caps, 12 or 8/strip	N801-0534 N801-0535
DNA Thermal Cycler 480 <sup>a</sup>	GeneAmp® Thin-Walled Reaction Tubes, 0.5-mL	N801-0537
	GeneAmp® Thin-Walled Reaction Tubes with Flat Cap	N801-0737
DNA Thermal Cycler (TC1) <sup>a</sup>	GeneAmp® Thin-Walled Reaction Tubes, 0.5-mL	N801-0537

a. The DNA Thermal Cycler (TC1) and the DNA Thermal Cycler 480 thermal cyclers require mineral oil that can be obtained from Applied Biosystems (P/N 0186-2302)



---

---

**Materials for  
Purifying  
Extension  
Products**

Method	Material	Supplier
Ethanol Precipitation	Ethanol (EtOH), non-denatured, 95%	MLS
	Aluminum foil tape, adhesive-backed	3M (Scotch Tape P/N 431 or 439) <sup>a</sup>

a. Contact 3M in the USA at (800) 364-3577 for your local 3M representative. Use of other tapes may result in leakage or contamination of the sample.

---

---

## Safety

---

---

**Documentation  
User Attention  
Words**

Five user attention words appear in the text of all Applied Biosystems user documentation. Each word implies a particular level of observation or action as described below.

**Note** Calls attention to useful information.

**IMPORTANT** Indicates information that is necessary for proper instrument operation.

**⚠ CAUTION** Indicates a potentially hazardous situation which, 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 which, if not avoided, could result in death or serious injury.

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

---

---

**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.

- 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.
- Minimize contact with and inhalation of chemicals. Wear appropriate personal protective equipment when handling

chemicals (e.g., safety glasses, gloves, or protective clothing). For additional safety guidelines, consult the MSDS.

- Do not leave chemical containers open. Use only with adequate ventilation.
- 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 Hazard Warning**

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

- 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.
- Handle chemical wastes in a fume hood.
- Minimize contact with and inhalation of chemical waste. Wear appropriate personal protective equipment when handling chemicals (e.g., safety glasses, gloves, or protective clothing).
- 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.

---

---

### **Site Preparation and Safety Guide**

A site preparation and safety guide is a separate document sent to all customers who have purchased an Applied Biosystems instrument. Refer to the guide written for your instrument for information on site preparation, instrument safety, chemical safety, and waste profiles.

---

---

### **About MSDSs**

Some of the chemicals used with this instrument may be listed as hazardous by their manufacturer. When hazards exist, warnings are prominently displayed on the labels of all chemicals.

Chemical manufacturers supply a current material safety data sheet (MSDS) before or with shipments of hazardous chemicals to new customers and with the first shipment of a hazardous chemical after an MSDS update. MSDSs provide you with the safety information you need to store, handle, transport and dispose of the chemicals safely.

We strongly recommend that you replace the appropriate MSDS in your files each time you receive a new MSDS packaged with a hazardous chemical.

**⚠ WARNING CHEMICAL HAZARD.** Be sure to familiarize yourself with the MSDSs before using reagents or solvents.

**Ordering MSDSs** You can order free additional copies of MSDSs for chemicals manufactured or distributed by Applied Biosystems using the contact information below.

To order MSDSs...	Then...							
Over the Internet	<p>a. Go to our Web site at <a href="http://www.appliedbiosystems.com/techsupp">www.appliedbiosystems.com/techsupp</a></p> <p>b. Click <b>MSDSs</b></p> <table border="1"> <thead> <tr> <th>If you have...</th> <th>Then...</th> </tr> </thead> <tbody> <tr> <td>The MSDS document number or the Document on Demand index number</td> <td>Enter one of these numbers in the appropriate field on this page.</td> </tr> <tr> <td>The product part number</td> <td rowspan="2">Select <b>Click Here</b>, then enter the part number or keyword(s) in the field on this page.</td> </tr> <tr> <td>Keyword(s)</td> </tr> </tbody> </table> <p>c. You can open and download a PDF (using Adobe® Acrobat® Reader™) of the document by selecting it, or you can choose to have the document sent to you by fax or e-mail.</p>	If you have...	Then...	The MSDS document number or the Document on Demand index number	Enter one of these numbers in the appropriate field on this page.	The product part number	Select <b>Click Here</b> , then enter the part number or keyword(s) in the field on this page.	Keyword(s)
If you have...	Then...							
The MSDS document number or the Document on Demand index number	Enter one of these numbers in the appropriate field on this page.							
The product part number	Select <b>Click Here</b> , then enter the part number or keyword(s) in the field on this page.							
Keyword(s)								
By automated telephone service	Use "To Obtain Documents on Demand" under "Technical Support."							
By telephone in the United States	Dial <b>1-800-327-3002</b> , then press <b>1</b> .							

To order MSDSs...	Then...	
By telephone from Canada	<b>To order in...</b>	<b>Dial 1-800-668-6913 and...</b>
	English	Press 1, then 2, then 1 again
	French	Press 2, then 2, then 1
By telephone from any other country	See the specific region under "To Contact Technical Support by Telephone or Fax" under "Technical Support."	

For chemicals not manufactured or distributed by Applied Biosystems, call the chemical manufacturer.

---

# *Preparing the Templates*

# 2

## Chapter Summary

---

**In This Chapter** The following topics are covered in this chapter:

<b>Topic</b>	<b>See Page</b>
Control DNA Templates	2-2
Template Preparation Methods	2-3
Single- and Double-Stranded Templates	2-3
BAC DNA Templates	2-3
PCR Templates	2-4
DNA Quantity	2-6

---

## Control DNA Templates

---

**Using Control DNA** Include a control DNA template as one of the templates in a set of sequencing reactions. The results from the control can help determine whether failed reactions are the result of poor template quality or sequencing reaction failure.

---

**Control DNA Sequence** We recommend M13mp18 as a single-stranded control and pGEM®-3Zf(+) as a double-stranded control. All Applied Biosystems DNA sequencing kits provide pGEM® control DNA.

The partial sequence of pGEM-3Zf(+) from the –21 M13 forward primer and M13 reverse primer, followed by the ensuing 1000 bases is shown in Appendix A, “Control DNA Sequences.”

---

**An Additional Control Sold Separately** The BigDye™ terminator v3.0 sequencing standard provides an additional control to help in troubleshooting electrophoresis runs. It contains lyophilized sequencing reactions that require only resuspension and denaturation before use.

There are two existing forms of the v3.0 sequencing standard, as shown in the table below. Please use the correct sequencing standard for your instrument. Refer to the product inserts for instructions on using each sequencing standard.

Instrument	Kit	Part Number
ABI PRISM® 3700 DNA Analyzer	ABI PRISM® 3700 BigDye™ Terminator v3.0 Sequencing Standard	4390309
ABI PRISM® 3100 Genetic Analyzer	ABI PRISM® BigDye™ Terminator v3.0 Sequencing Standard	4390303
ABI PRISM® 310 Genetic Analyzer		
ABI PRISM® 377 DNA Sequencers <sup>a</sup>		
ABI PRISM® 373 DNA Sequencers with the BigDye™ Filter Wheel <sup>b</sup>		

a. Includes the ABI PRISM 377, ABI PRISM 377-18, ABI PRISM 377 with XL Upgrade, and the ABI PRISM 377 with 96-Lane Upgrade instruments.

b. Includes the ABI PRISM 373 and ABI PRISM 373 with XL Upgrade instruments.

## Template Preparation Methods

---

**Single- and Double-Stranded Templates** Refer to the *Automated DNA Sequencing Chemistry Guide* (P/N 4305080) for information on preparing single- and double-stranded templates.

---

**BAC DNA Templates** With larger DNA targets such as bacterial artificial chromosomes (BACs), the quality of DNA template is important to the success of the sequencing reaction. Two methods have given good sequencing results:

- Alkaline lysis<sup>1</sup>, with extra phenol extraction and isopropanol precipitation if very clean DNA is desired
- Cesium chloride (CsCl) banding

### Commercial Kits

Commercial kits are also available for BAC DNA preparation:

- QIAGEN-tip 100 (QIAGEN: P/N 10043, 25 reactions; 10045, 100 reactions)
  - QIAGEN-tip 500 (QIAGEN: P/N 10063, 25 reactions; 10065, 100 reactions)
- 

1. Marra, M., Weinstock, L.A., and Mardis, E.R. 1996. End sequence determination from large insert cloning using energy transfer fluorescent primers. *Genomic Methods* 6: 1118–1122.

---

---

**PCR Templates** Cycle sequencing provides the most reproducible results for sequencing PCR templates. Although PCR fragments can be difficult to denature with traditional sequencing methods, cycle sequencing provides several chances to denature and extend the template, which ensures adequate signal in the sequencing reaction.

**Importance of Purifying Product**

For optimum results, purify the PCR product before sequencing. In general, any method that removes dNTPs and primers should work. We recommend Centricon®-100 columns (P/N N930-2119). The protocol for using these columns is provided in “Purifying PCR Fragments” on page 2-5.

Refer to the *Automated DNA Sequencing Chemistry Guide* (P/N 4305080) for information on sequencing PCR templates.

---

---

**Direct Link PCR Sequencing** PCR protocols that limit amounts of primers and dNTPs allow the product of the reaction to be used for sequencing with no purification. This is usually carried out by setting up the PCR amplification with 5–10 pmol of primers and 20–40 μM dNTPs, so that most of the primers and dNTPs are exhausted during amplification.

If the yield of the desired PCR product is high and the product is specific, *i.e.*, it produces a single band when analyzed by agarose gel electrophoresis, the sample can be diluted before sequencing and will give good results. The dilution ratio depends on the concentration of your PCR product and needs to be determined empirically (start with 1:2 and 1:10 dilutions with deionized water).

When you limit concentrations of primers and dNTPs and dilute the PCR products, the PCR parameters have to be robust. Direct Link PCR sequencing is most useful in applications where the same target is being amplified and sequenced repeatedly and PCR conditions have been optimized.

---

---



---

---

**Purifying PCR  
Fragments**

To purify PCR fragments by ultrafiltration:

<b>Step</b>	<b>Action</b>
<b>1</b>	Assemble the Centricon-100 column according to the manufacturer's recommendations.
<b>2</b>	Load 2 mL deionized water onto the column.
<b>3</b>	Add the entire sample to the column.
<b>4</b>	Spin the column at $3000 \times g$ in a fixed-angle centrifuge for 10 minutes.  <b>Note</b> The manufacturer recommends a maximum speed of $1000 \times g$ , but $3000 \times g$ has worked well in Applied Biosystems laboratories. If you are following the manufacturer's guidelines, increase the time to compensate.
<b>5</b>	Remove the waste receptacle and attach the collection vial.
<b>6</b>	Invert the column and spin it at $270 \times g$ for 2 minutes to collect the sample. This should yield approximately 40–60 $\mu\text{L}$ of sample.
<b>7</b>	Add deionized water to bring the purified PCR fragments to the original volume.

---

---

## DNA Quantity

---

**Quantitating DNA** For most purified PCR products, using 1  $\mu\text{L}$  for each reaction (A, C, G, and T) will work. If possible, quantitate the amount of purified DNA by measuring the absorbance at 260 nm or by some other method. If you do not quantitate the amount of DNA beforehand, you run the risk of using too much DNA, which can affect the final sequencing result.

---

**Template Quantity** The table below shows the amount of template to use in a cycle sequencing reaction.

Template	Quantity
PCR product:	
100–200 bp	2–5 ng
200–500 bp	5–10 ng
500–1000 bp	10–20 ng
1000–2000 bp	20–50 ng
>2000 bp	50–150 ng
Single-stranded	100–400 ng
Double-stranded	200–800 ng
Cosmid, BAC	0.5–1.0 $\mu\text{g}$

**Note** In general, higher DNA quantities give higher signal intensities.

The ranges given in the table above should work for all primers. You may be able to use even less DNA, especially when sequencing with the –21 M13 primer. The amount of PCR product to use in sequencing will also depend on the length and purity of the PCR product.

---

# *Performing Cycle Sequencing*

# 3

## Chapter Summary

---

**In This Chapter** The following topics are covered in this chapter:

<b>Topic</b>	<b>See Page</b>
Introduction	3-2
Cycle Sequencing Single- and Double-Stranded DNA	3-3
Cycle Sequencing BAC DNA	3-8

---

## Introduction

---

**Overview** The cycle sequencing protocols used for the ABI PRISM® BigDye™ Primer v3.0 Ready Reaction Cycle Sequencing Kit with AmpliTaq® DNA Polymerase, FS, are the same as those used for the ABI PRISM® BigDye™ Primer original Ready Reaction Cycle Sequencing Kit. They have been optimized for Applied Biosystems thermal cyclers.

---

**Optimizing Cycle Parameters** The standard cycling conditions used in this manual have been optimized for the –21 M13 forward primer and work well for most primers. If you obtain low signal with the M13 reverse primer, you can improve your sequencing data by altering the cycling conditions as follows:

- Standard conditions use 55 °C as the annealing temperature in the three-step cycle. Lower the annealing temperature to 45–48 °C.
  - The number of three-step versus two-step cycles can also affect signal strength. Increase the number of three-step cycles to 20 instead of the standard 15 (increasing the total number of cycles to 35).
-

## Cycle Sequencing Single- and Double-Stranded DNA

---

**Overview** This section describes how to prepare reactions and perform cycle sequencing on a variety of templates, including M13, plasmids, and PCR products.

---

**Preparing the Reactions** **Three Options**  
The flexibility of the BigDye™ primers v3.0 allows three options for cycle sequencing:

Reaction Type	Template	Cycle
1X	<ul style="list-style-type: none"><li>• PCR product</li><li>• plasmid</li><li>• M13</li></ul>	standard
0.5X	<ul style="list-style-type: none"><li>• PCR product</li><li>• plasmid</li><li>• M13</li></ul>	standard
High-sensitivity (2X)	<ul style="list-style-type: none"><li>• BACs</li><li>• extra long PCR products</li><li>• other large DNA</li></ul>	modified

### Procedures

The type of tube required depends on the thermal cycler that you are using. Refer to “Materials for Cycle Sequencing” on page 1-10.

**IMPORTANT** Prepare separate tubes for each of the four reactions (A, C, G, and T).

### 1X Reactions

Step	Action																				
1	Aliquot the following reagents into four PCR tubes:																				
	<table border="1"> <thead> <tr> <th>Reagent</th> <th>A (μL)</th> <th>C (μL)</th> <th>G (μL)</th> <th>T (μL)</th> </tr> </thead> <tbody> <tr> <td>Ready Reaction Premix</td> <td>4</td> <td>4</td> <td>4</td> <td>4</td> </tr> <tr> <td>DNA Template (see page 2-6 for quantity)</td> <td>1</td> <td>1</td> <td>1</td> <td>1</td> </tr> <tr> <td>Total Volume</td> <td>5</td> <td>5</td> <td>5</td> <td>5</td> </tr> </tbody> </table>	Reagent	A (μL)	C (μL)	G (μL)	T (μL)	Ready Reaction Premix	4	4	4	4	DNA Template (see page 2-6 for quantity)	1	1	1	1	Total Volume	5	5	5	5
	Reagent	A (μL)	C (μL)	G (μL)	T (μL)																
	Ready Reaction Premix	4	4	4	4																
DNA Template (see page 2-6 for quantity)	1	1	1	1																	
Total Volume	5	5	5	5																	
2	<b>If using the DNA Thermal Cycler (TC1) or DNA Thermal Cycler 480:</b>																				
	a. Add 20 μL of light mineral oil. b. Spin to layer the oil over the aqueous reaction.																				

### 0.5X Reactions

Dilute 5X Sequencing Buffer (P/N 4305605, 600 reactions; 4305603, 5400 reactions) with four parts deionized water to 1X for use in this procedure.

Step	Action																				
1	Dilute each Ready Reaction Premix (A, C, G, T) 1:1 with 1X Sequencing Buffer in a separate tube (e.g., 2 μL of A Mix and 2 μL of 1X Sequencing Buffer).																				
2	Aliquot the following reagents into four PCR tubes for each DNA template:																				
	<table border="1"> <thead> <tr> <th>Reagent</th> <th>A (μL)</th> <th>C (μL)</th> <th>G (μL)</th> <th>T (μL)</th> </tr> </thead> <tbody> <tr> <td>Diluted Ready Reaction Premix</td> <td>4</td> <td>4</td> <td>4</td> <td>4</td> </tr> <tr> <td>DNA Template (see page 2-6 for quantity)</td> <td>1</td> <td>1</td> <td>1</td> <td>1</td> </tr> <tr> <td>Total Volume</td> <td>5</td> <td>5</td> <td>5</td> <td>5</td> </tr> </tbody> </table>	Reagent	A (μL)	C (μL)	G (μL)	T (μL)	Diluted Ready Reaction Premix	4	4	4	4	DNA Template (see page 2-6 for quantity)	1	1	1	1	Total Volume	5	5	5	5
	Reagent	A (μL)	C (μL)	G (μL)	T (μL)																
	Diluted Ready Reaction Premix	4	4	4	4																
DNA Template (see page 2-6 for quantity)	1	1	1	1																	
Total Volume	5	5	5	5																	
3	<b>If using the DNA Thermal Cycler (TC1) or DNA Thermal Cycler 480:</b>																				
	a. Add 20 μL of light mineral oil. b. Spin to layer the oil over the aqueous reaction.																				

### High-Sensitivity (2X) Reactions

Aliquot the following reagents into four PCR tubes:

Reagent	A (µL)	C (µL)	G (µL)	T (µL)
Ready Reaction Premix	8	8	8	8
DNA Template (see page 2-6 for quantity)	2	2	2	2
Total Volume	10	10	10	10

**Note** Use the GeneAmp® PCR System 9600 or 9700 for cycle sequencing high-sensitivity reactions. The use of other thermal cyclers is not supported.

If the majority of your applications require high sensitivity, please contact your local Field Applications Specialist or Technical Support for more options.

---

---

---

**Cycle Sequencing  
on the System  
9700, 9600, or 2400**

**Note** The total time required for the cycling in this protocol is approximately 1 hour and 15 minutes.

To sequence single- and double-stranded DNA on the GeneAmp® PCR System 9700 (in 9600 emulation mode), 9600, or 2400:

<b>Step</b>	<b>Action</b>
<b>1</b>	Place the tubes in a thermal cycler and set the volume to 5 µL.
<b>2</b>	Repeat the following for 15 cycles: <ul style="list-style-type: none"><li>• Rapid thermal ramp<sup>a</sup> to 96 °C</li><li>• 96 °C for 10 sec.</li><li>• Rapid thermal ramp to 55 °C</li><li>• 55 °C for 5 sec.</li><li>• Rapid thermal ramp to 70 °C</li><li>• 70 °C for 1 min.</li></ul>
<b>3</b>	Repeat the following for 15 cycles: <ul style="list-style-type: none"><li>• Rapid thermal ramp to 96 °C</li><li>• 96 °C for 10 sec.</li><li>• Rapid thermal ramp to 70 °C</li><li>• 70 °C for 1 min.</li></ul>
<b>4</b>	Rapid thermal ramp to 4 °C and hold until ready to precipitate.

a. Rapid thermal ramp is 1 °C/sec.

---

---



---

---

**Cycle Sequencing  
on the TC1 or 480**

**Note** The total time required for the cycling in this protocol is approximately 1 hour and 45 minutes.

To sequence single- and double-stranded on the DNA Thermal Cycler (TC1) or DNA Thermal Cycler 480:

<b>Step</b>	<b>Action</b>
<b>1</b>	Place the tubes in a thermal cycler preheated to 95 °C.
<b>2</b>	Repeat the following for 15 cycles: <ul style="list-style-type: none"><li>• Rapid thermal ramp<sup>a</sup> to 95 °C</li><li>• 95 °C for 30 sec.</li><li>• Rapid thermal ramp to 55 °C</li><li>• 55 °C for 30 sec.</li><li>• Rapid thermal ramp to 70 °C</li><li>• 70 °C for 1 min.</li></ul>
<b>3</b>	Repeat the following for 15 cycles: <ul style="list-style-type: none"><li>• Rapid thermal ramp to 95 °C</li><li>• 95 °C for 30 sec.</li><li>• Rapid thermal ramp to 70 °C</li><li>• 70 °C for 1 min.</li></ul>
<b>4</b>	Rapid thermal ramp to 4 °C and hold until ready to precipitate.

a. Rapid thermal ramp is 1 °C/sec.

---

---

## Cycle Sequencing BAC DNA

---

**Preparing Sequencing Reactions** Follow the procedures given for preparing reactions in “Cycle Sequencing Single- and Double-Stranded DNA” (see “Preparing the Reactions” on page 3-3).

---

**Performing Cycle Sequencing** **Note** The total time required for the cycling in this protocol is approximately 1 hour and 45 minutes.

To sequence BAC DNA on the GeneAmp PCR System 9700 (in 9600 emulsion mode), 9600, or 2400:

Step	Action
1	Place the tubes in a thermal cycler, set the volume at 10 $\mu$ L, and begin thermal cycling with the following parameters: <ul style="list-style-type: none"><li>• Rapid thermal ramp<sup>a</sup> to 95 °C</li><li>• 95 °C for 5 min.</li></ul>
2	Repeat the following for 20 cycles: <ul style="list-style-type: none"><li>• Rapid thermal ramp to 95 °C</li><li>• 95 °C for 30 sec.</li><li>• Rapid thermal ramp to 50 °C</li><li>• 50 °C for 15 sec.</li><li>• Rapid thermal ramp to 70 °C</li><li>• 70 °C for 1 min.</li></ul>
3	Repeat the following for 15 cycles: <ul style="list-style-type: none"><li>• Rapid thermal ramp to 95 °C</li><li>• 95 °C for 30 sec.</li><li>• Rapid thermal ramp to 70 °C</li><li>• 70 °C for 1 min.</li></ul>
4	Rapid thermal ramp to 4 °C and hold until ready to precipitate.

a. Rapid thermal ramp is 1 °C/sec.

---

# *Preparing the Samples for Loading*

# 4

## Chapter Summary

---

**In This Chapter** The following topics are covered in this chapter:

Topic	See Page
Recommendations for Precipitation/Loading	4-2
Ethanol Precipitation for Microcentrifuge Tubes	4-3
Ethanol Precipitation for 96-Well Reaction Plates	4-5

---

## Recommendations for Precipitation/Loading

---

**For the 3700, 3100, 310, and 373 Instruments** In the ethanol precipitation method, reactions from four separate tubes are pooled together and precipitated using ethanol.

This method is recommended for the following instruments:

- ABI PRISM® 3700 DNA Analyzer
- ABI PRISM® 3100 Genetic Analyzer
- ABI PRISM® 310 Genetic Analyzer
- ABI PRISM® 373 DNA Sequencers with the ABI PRISM® BigDye™ Filter Wheel installed.<sup>1</sup>

For the 3700, 3100, and 310 instruments, use this method to remove salts and other components of the reaction. For the 373 instruments, use this method to concentrate the samples.

See “Ethanol Precipitation for Microcentrifuge Tubes” on page 4-3 or “Ethanol Precipitation for 96-Well Reaction Plates” on page 4-5.

---

**For the 377 Instrument** For the ABI PRISM® 377 DNA Sequencer (all models<sup>2</sup>), there are two options:

- Ethanol precipitation—samples are concentrated.  
See “Ethanol Precipitation for Microcentrifuge Tubes” on page 4-3 or “Ethanol Precipitation for 96-Well Reaction Plates” on page 4-5.
- Express Load—samples are loaded directly onto a gel without being precipitated  
See “Express Load Option for 36-Well Gels” on page 5-11.

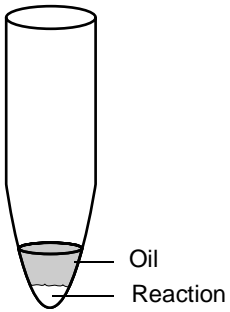
**IMPORTANT** The Express Load option is only for the 377 instrument. It cannot be used for the 3700, 3100, 310, or 373 instruments.

---

1. Includes the ABI PRISM 373 and ABI PRISM 373 with XL Upgrade instruments.
2. Includes the ABI PRISM 377, ABI PRISM 377-18, ABI PRISM 377 with XL Upgrade, and the ABI PRISM 377 with 96-Lane Upgrade instruments.

## Ethanol Precipitation for Microcentrifuge Tubes

**Procedure** To precipitate in microcentrifuge tubes:

Step	Action
1	<p>To a clean microcentrifuge tube, add:</p> <ul style="list-style-type: none"> <li>• 60 <math>\mu</math>L of non-denatured 95% ethanol</li> </ul> <p>OR</p> <ul style="list-style-type: none"> <li>• 120 <math>\mu</math>L of non-denatured 95% ethanol, if sequencing BAC DNA or other high-sensitivity reactions</li> </ul> <p><b>Note</b> The use of sodium acetate is not necessary for precipitation.</p> <p><b>⚠ WARNING CHEMICAL HAZARD.</b> Ethanol is a flammable liquid and vapor. It may cause eye, skin, and upper respiratory tract irritation. Prolonged or repeated contact may dry skin. Exposure may cause central nervous system depression and liver damage. Please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.</p>
2	<p>Pipet the extension reactions from the bottom of each of the four tubes into the ethanol mixture. Mix thoroughly.</p> <p><b>IMPORTANT To remove reactions run on the TC1 or DNA Thermal Cycler 480:</b> Place the pipette tip into the bottom of the reaction and carefully remove the reaction from the oil (see “Procedural Options” on page 4-4). Transfer as little oil as possible.</p>  <p>The diagram shows a microcentrifuge tube with a small amount of liquid at the bottom. A horizontal line indicates the interface between two layers. The upper layer is labeled 'Oil' and the lower layer is labeled 'Reaction'.</p>
3	<p>Place the tube on wet ice or leave it at room temperature for 10–15 minutes to precipitate the extension products.</p>
4	<p>Spin the tube in a microcentrifuge for 10–20 minutes at maximum speed. Carefully aspirate or decant the supernatant and discard. At this point, a pellet may or may not be visible.</p>

To precipitate in microcentrifuge tubes: *(continued)*

Step	Action
5	<p><b>Optional:</b> Rinse the pellet with 250 <math>\mu\text{L}</math> of 70% ethanol and spin for 5 minutes in a microcentrifuge. Again, carefully aspirate or decant the supernatant and discard. This may remove some of the salts from the pellet, but doing so is often not necessary.</p> <p><b>Note</b> If you use sodium acetate, you must rinse the pellet. This will reduce the carryover of salt.</p>
6	Dry the pellet in a vacuum centrifuge for 1–3 minutes, or until dry. Do not overdry.

---

**Procedural Options** Removing the four reactions from under the oil can be difficult because of their small (5  $\mu\text{L}$ ) volumes. If this is a problem, try one of the following options:

- After cycling, add 5  $\mu\text{L}$  of deionized water to each reaction for a total volume of 10  $\mu\text{L}$ .  
If this is done, increase the amount of 95% ethanol used in step 1 of “Ethanol Precipitation for Microcentrifuge Tubes” on page 4-3 to 120  $\mu\text{L}$ .
  - Double the amount of Ready Reaction Premix and DNA template for each reaction (A, C, G, and T) to yield several loadings’ worth (*i.e.*, a 2X reaction).
-

## Ethanol Precipitation for 96-Well Reaction Plates

**Standard Method** With the BigDye™ primers v3.0, the standard method is ethanol precipitation without the addition of salt. A final 70% ethanol wash is optional.

**IMPORTANT** Use non-denatured 95% ethanol rather than absolute (100%) ethanol. Absolute ethanol absorbs water from the atmosphere, gradually decreasing its concentration. This can lead to inaccurate final concentrations of ethanol, which can affect some protocols.

### Precipitating in 96-Well Reaction Plates

To precipitate in 96-well reaction plates:

Step	Action
1	Remove the four 96-well reaction plates from the thermal cycler. Remove the caps from each tube.
2	Briefly centrifuge the plates to collect the samples at the bottoms of the tubes.
3	Using a multichannel pipet, pool the entire samples' volumes from each of the four plates ( <i>i.e.</i> , containing the dideoxy A, C, G, and T reactions).
4	To the pooled reactions, add: <ul style="list-style-type: none"><li>• 60 µL of non-denatured 95% ethanol (EtOH)</li></ul> <p><b>⚠ WARNING CHEMICAL HAZARD.</b> Ethanol is a flammable liquid and vapor. It may cause eye, skin, and upper respiratory tract irritation. Prolonged or repeated contact may dry skin. Exposure may cause central nervous system depression and liver damage. Please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.</p>
5	Seal the tubes with strip caps or by applying a piece of 3M Scotch Tape 431 or 439 adhesive-backed aluminum foil tape. Press the foil onto the tubes to prevent any leakage.
6	Invert the plate a few times to mix.
7	Leave the plate on ice or at 4 °C for 15 minutes to precipitate the extension products.

To precipitate in 96-well reaction plates: *(continued)*

Step	Action
8	<p>Place the plate in a table-top centrifuge with a plate adaptor and spin it at the maximum speed, which must be <math>\geq 1400 \times g</math> but <math>&lt; 3000 \times g</math>:</p> <ul style="list-style-type: none"><li>• <math>1400\text{--}2000 \times g</math>: 45 minutes</li><li>• <math>2000\text{--}3000 \times g</math>: 30 minutes</li></ul> <p><b>IMPORTANT</b> Proceed to the next step immediately. If this is not possible, then spin the tubes for 2 minutes more immediately before performing the next step.</p>
9	<p>Without disturbing the precipitates, remove the adhesive tape and discard the supernatant by inverting the plate onto a paper towel folded to the size of the plate.</p>
10	<p>Place the inverted plate with the paper towel into the table-top centrifuge and spin at <math>50 \times g</math> for 1 minute.</p>
11	<p>Remove the plate and discard the paper towel.</p> <p><b>Note</b> Pellets may or may not be visible. Vacuum drying of the samples is not necessary.</p>

---

---



# *Sample Electrophoresis*

# 5

## Chapter Summary

---

**In This Chapter** The following topics are covered in this chapter:

<b>Topic</b>	<b>See Page</b>
Before You Begin	5-2
Electrophoresis on the ABI PRISM 3700 DNA Analyzer	5-3
Electrophoresis on the ABI PRISM 3100 Genetic Analyzer	5-4
Electrophoresis on the ABI PRISM 310 Genetic Analyzer	5-5
Electrophoresis on the ABI PRISM 377 DNA Sequencers	5-8
Electrophoresis on the 373 Instrument with BigDye Filter Wheel	5-8

---

## Before You Begin

- 
- Important Reminders**
- Be sure to use the correct dye set/primer (mobility) file for the BigDye™ primers v3.0. The dye base relationships are different than with the BigDye™ terminators v3.0 and dRhodamine terminators.
  - For the ABI PRISM® 3700 DNA Analyzer, ABI PRISM® 3100 Genetic Analyzer, and ABI PRISM® 310 Genetic Analyzer, separate dye set/primer (mobility) files are used for the forward and reverse primers.
  - We do not recommend using POP-5™ polymer on the 3700 instrument for BigDye primer v3.0 chemistry.
  - For the ABI PRISM® 377 DNA Sequencers, a single dye set/primer (mobility) file is used for the forward and reverse primers.
-

## Electrophoresis on the ABI PRISM 3700 DNA Analyzer

---

**Requirements** Electrophoresis and data analysis of samples on the ABI PRISM® 3700 DNA Analyzer requires the following:

### Run Modules

Configuration	Run Module
POP-6™ polymer, 50-cm	Seq1_1POP6DefaultModule

### Dye Set/Primer (Mobility) Files

Polymer	Dye Set/Primer (Mobility) File
POP-6 polymer	DP3700POP6{BDv3-21M13}v1.mob
	DP3700POP6{BDv3-21M13Rev}v1.mob

**Note** For the 3700 instrument, separate dye set/primer (mobility) files are used for the forward and reverse primers.

### Standards

**IMPORTANT** Use Dye Set D.

Dye Set	Standards for Spectral Calibration
D	ABI PRISM® 3700 BigDye™ Terminator v3.0 Sequencing Standard (P/N 4390309)

**Note** Refer to the product insert for instructions on using the standards for this instrument.

---

**Performing Sample Electrophoresis** For information on how to perform sample electrophoresis on the 3700 instrument, refer to the following manuals:

- *ABI PRISM 3700 DNA Analyzer Sequencing Chemistry Guide* (P/N 4309125)
- *ABI PRISM 3700 DNA Analyzer User's Manual* (P/N 4306152)

---

## Electrophoresis on the ABI PRISM 3100 Genetic Analyzer

---

**Requirements** Electrophoresis and data analysis of samples on the ABI PRISM® 3100 Genetic Analyzer requires the following:

### Run Modules

Configuration	Run Module
POP-6™ polymer, 36-cm	RapidSeq36_POP6DefaultModule
POP-6 polymer, 50-cm	StdSeq50_POP6DefaultModule

### Dye Set/Primer (Mobility) Files

Polymer	Dye Set/Primer (Mobility) File
POP-6 polymer	DP3100POP6{BDv3-21M13}v1.mob
	DP3100POP6{BDv3-21M13Rev}v1.mob

**Note** For the 3100 instrument, separate dye set/primer (mobility) files are used for the forward and reverse primers.

### Standards

**IMPORTANT** Use Dye Set Z.

Dye Set	Standards for Spectral Calibration
Z	ABI PRISM® BigDye™ Terminator v3.0 Sequencing Standard (P/N 4390303)

**Note** Refer to the product insert for instructions on using the standards for this instrument.

---

**Performing Sample Electrophoresis** For information on how to perform sample electrophoresis on the 3100 instrument, refer to the following manuals:

- *ABI PRISM 3100 Genetic Analyzer Sequencing Chemistry Guide* (P/N 4315831)
- *ABI PRISM 3100 Genetic Analyzer User's Manual* (P/N 4315834)

---

## Electrophoresis on the ABI PRISM 310 Genetic Analyzer

**Requirements** Electrophoresis and data analysis of samples on the ABI PRISM® 310 Genetic Analyzer requires the following:

### Filter Set E Run Modules

Configuration	Run Module
POP-4™ polymer, 1-mL syringe, 47-cm, 50-µm i.d. capillary, Ld = 36 cm	P4StdSeq (1 mL) E
POP-4 polymer, Rapid Sequencing, 1-mL syringe, 47-cm, 50-µm i.d. capillary, Ld = 36 cm	P4RapidSeq (1 mL) E
POP-6™ polymer, 1-mL syringe, 61-cm, 50-µm i.d. capillary	Seq POP6 (1 mL) E
POP-6 polymer, Rapid Sequencing, 1-mL syringe, 47-cm, 50-µm i.d. capillary	Seq POP6 Rapid (1 mL) E

### Dye Set/Primer (Mobility) Files

Polymer	Dye Set/Primer (Mobility) File
POP-4 polymer	DP310POP4{BDv3-21M13}v1.mob
	DP310POP4{BDv3-M13Rev}v1.mob
POP-6 polymer	DP310POP6{BDv3-21M13}v1.mob
	DP310POP6{BDv3-M13Rev}v1.mob

**Note** For the 310 instrument, separate dye set/primer (mobility) files are used for the forward and reverse primers.

### Matrix Standards

**IMPORTANT** The instrument (matrix) file for the BigDye primers v3.0 cannot be used for the BigDye terminators (original), BigDye terminators v2.0, dRhodamine terminators, or BigDye primers (original).

Dye/Filter Set	Standards for Instrument (Matrix) File Generation
E	ABI PRISM® BigDye™ Matrix Standards v3.0 (P/N 4390421)

**Note** Refer to the product insert for instructions on using the standards for this instrument.

---

**Setting the Data Analysis Range** The unincorporated dye primer peak at the beginning of the electropherogram (and the full-length PCR peak at the end of the electropherogram if sequencing PCR products) must be excluded from analysis for data to be normalized correctly.

Analysis of ABI PRISM 310 instrument data is especially sensitive to the location of the Peak 1 Location and Analysis Start Point. The software should pick these correctly in most cases. A miscalled Peak 1 Location and/or Analysis Start Point can result in spacing anomalies visible in the analyzed data.

Refer to page 6-15 of the *Automated DNA Sequencing Chemistry Guide* (P/N 4305080) for more information on setting the data analysis range.

---

**Resuspending the Samples for Sequencing with POP-6 Polymer** To resuspend the samples for sequencing with POP-6 polymer:

Step	Action
1	Resuspend each sample pellet in 25 $\mu$ L of template suppression reagent (TSR, supplied with the polymer).  <b>⚠ CAUTION CHEMICAL HAZARD. Template suppression reagent</b> may cause eye, skin, and respiratory tract irritation. Please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.
2	Vortex and spin the samples.
3	Heat the samples at 95 °C for 2 minutes to denature, then chill on ice.
4	Vortex and spin the samples again. Place on ice until ready to use.
5	Refer to the <i>ABI PRISM 310 Genetic Analyzer User's Manual</i> (P/N 903565) for guidelines on loading the samples.

**Note** Although not recommended on a routine basis, you can keep samples prepared in TSR frozen for several weeks before running on the ABI PRISM 310 Genetic Analyzer with no detectable loss in resolution or base calling.

---

---

---

**Resuspending the  
Samples for  
Sequencing with  
POP-4 Polymer**

Deionized formamide is the recommended resuspension solution for sequencing with POP-4 polymer on the 310 genetic analyzer.

**IMPORTANT** Use only the highest grade of deionized formamide, such as Applied Biosystems Hi-Di™ formamide (P/N 4311320). Deionized formamide should be stored at –20 °C in usable aliquots to prevent several freeze-thaw cycles. If the deionized formamide stored at –20 °C does not freeze, discard it and use fresh deionized formamide for sample resuspension.

To resuspend the samples for sequencing with POP-4 polymer:

Step	Action
1	Resuspend each sample pellet in 25 µL of deionized formamide.  <b>IMPORTANT</b> Securely seal each sample tube after resuspension with deionized formamide to limit the sample's exposure to air.  <b>⚠ WARNING CHEMICAL HAZARD. Formamide</b> is harmful if absorbed through the skin and may cause irritation to the eyes, skin, and respiratory tract. It may cause damage to the central nervous system and the male and female reproductive systems, and is a possible birth defect hazard. Please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.
2	Vortex and spin the samples.
3	Heat the samples at 95 °C for 2 minutes to denature, then chill on ice.
4	Vortex and spin the samples again. Place on ice until ready to use.
5	Refer to the <i>ABI PRISM 310 Genetic Analyzer User's Manual</i> (P/N 903565) for guidelines on loading the samples.

**Note** Although extended sample storage at room temperature is not recommended on a routine basis, you can keep securely sealed samples prepared in deionized formamide at room temperature for up to 48 hours on the ABI PRISM 310 Genetic Analyzer autosampler with no detectable loss in resolution or base calling. Freezing of samples resuspended in deionized formamide is not advised.

---

---

## Electrophoresis on the ABI PRISM 377 DNA Sequencers

**Requirements** Electrophoresis and data analysis of samples on the ABI PRISM® 377 DNA Sequencers (all models<sup>1</sup>) require the following:

### Filter Set E Run Modules

Configuration <sup>a</sup>	Run Module
36-cm wtr, 1200 scans/hr, any comb	Seq Run 36E-1200
36-cm wtr, 2400 scans/hr, any comb	Seq Run 36E-2400
48-cm wtr, 1200 scans/hr, any comb	Seq Run 48E-1200

a. Any plate check and prerun module can be used on the ABI PRISM 377 DNA Sequencers.

### Dye Set/Primer (Mobility) Files

Gel Formulation	Dye Set/Primer (Mobility) File
4.5% acrylamide (29:1) or 5% Long Ranger™ gel	DP377{BDv3-21M13}v1.mob
	DP377{BDv3-M13Rev}v1.mob

**Note** For the ABI PRISM 377 DNA Sequencers, a single dye set/primer (mobility) file is used for the forward and reverse primers.

### Matrix Standards

**IMPORTANT** The instrument (matrix) file for the BigDye primers v3.0 cannot be used for the BigDye terminators (original), BigDye terminators v2.0, dRhodamine terminators, or BigDye primers (original).

Dye/Filter Set	Standards for Instrument (Matrix) File Generation
E	ABI PRISM® BigDye™ Matrix Standards v3.0 (P/N 4390421)

**Note** Refer to the product insert for instructions on using the standards for this instrument.

---

1. Includes the ABI PRISM 377, ABI PRISM 377-18, ABI PRISM 377 with XL Upgrade, and the ABI PRISM 377 with 96-Lane Upgrade instruments.



---

---

**Using the Lane Guide Kit** If you are using the BigDye chemistries v3.0 on the 377 instrument in conjunction with the ABI PRISM® Lane Guide™ Lane Identification Kit, refer to that kit's protocol (P/N 4313804) for instructions on resuspending and loading samples.

---

---

**Using Long-Read Gel and Buffer Formulations** For longer sequencing read lengths follow the gel and buffer formulations described in the user bulletin entitled *Achieving Longer High Accuracy Reads on the 377 Sequencer* (P/N 4315153).

---

---

**Resuspending and Loading the Samples**

**Running Ethanol-Precipitated Samples**

Step	Action															
1	<p>Prepare a loading buffer by combining the following in a 5:1 ratio (5 parts deionized formamide to 1 part EDTA with blue dextran):</p> <ul style="list-style-type: none"> <li>• Deionized formamide</li> <li>• 25 mM EDTA (pH 8.0) with blue dextran (50 mg/mL)</li> </ul> <p><b>⚠ WARNING CHEMICAL HAZARD.</b> Formamide is harmful if absorbed through the skin and may cause irritation to the eyes, skin, and respiratory tract. It may cause damage to the central nervous system and the male and female reproductive systems, and is a possible birth defect hazard. Please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.</p> <p><b>⚠ CAUTION CHEMICAL HAZARD.</b> EDTA may cause eye, skin, and respiratory tract irritation. Please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.</p>															
2	<p>Resuspend each sample pellet in loading buffer as follows:</p> <table border="1"> <thead> <tr> <th>Template</th> <th>Reaction</th> <th>Volume (μL), 36-well</th> <th>Volume (μL), 48- or 64-well</th> </tr> </thead> <tbody> <tr> <td rowspan="2">PCR product, plasmid, M13</td> <td>1X</td> <td>6</td> <td>4</td> </tr> <tr> <td>0.5X</td> <td>2–4</td> <td>1–2</td> </tr> <tr> <td>BAC, large DNA</td> <td>2X</td> <td>2–4</td> <td>1–2</td> </tr> </tbody> </table>	Template	Reaction	Volume (μL), 36-well	Volume (μL), 48- or 64-well	PCR product, plasmid, M13	1X	6	4	0.5X	2–4	1–2	BAC, large DNA	2X	2–4	1–2
Template	Reaction	Volume (μL), 36-well	Volume (μL), 48- or 64-well													
PCR product, plasmid, M13	1X	6	4													
	0.5X	2–4	1–2													
BAC, large DNA	2X	2–4	1–2													
3	Vortex and spin the samples.															
4	Heat the samples at 95 °C for 2 minutes to denature. Place on ice until ready to load.															
5	<p>Load each sample into a separate lane of the gel as follows:</p> <table border="1"> <thead> <tr> <th>Template</th> <th>Reaction</th> <th>Volume (μL), 36-well</th> <th>Volume (μL), 48- or 64-well</th> </tr> </thead> <tbody> <tr> <td rowspan="2">PCR product, plasmid, M13</td> <td>1X</td> <td>0.75–2.0</td> <td>0.5–1.5</td> </tr> <tr> <td>0.5X</td> <td>2–3</td> <td>1–1.5</td> </tr> <tr> <td>BAC, large DNA</td> <td>2X</td> <td>2–3</td> <td>1–1.5</td> </tr> </tbody> </table>	Template	Reaction	Volume (μL), 36-well	Volume (μL), 48- or 64-well	PCR product, plasmid, M13	1X	0.75–2.0	0.5–1.5	0.5X	2–3	1–1.5	BAC, large DNA	2X	2–3	1–1.5
Template	Reaction	Volume (μL), 36-well	Volume (μL), 48- or 64-well													
PCR product, plasmid, M13	1X	0.75–2.0	0.5–1.5													
	0.5X	2–3	1–1.5													
BAC, large DNA	2X	2–3	1–1.5													

### Express Load Option for 36-Well Gels

Step	Action
1	<p>Combine the four reactions (A, C, G, T) with 5 <math>\mu\text{L}</math> of 5 mM EDTA (25 <math>\mu\text{L}</math> total volume).</p> <p><b>⚠ CAUTION CHEMICAL HAZARD.</b> EDTA may cause eye, skin, and respiratory tract irritation. Please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.</p>
2	<p>Vortex briefly, then spin in a microcentrifuge.</p>
3	<p>Prepare a loading buffer by combining the following in a 5:1 ratio (5 parts deionized formamide to 1 part EDTA with blue dextran):</p> <ul style="list-style-type: none"><li>• Deionized formamide</li><li>• 25 mM EDTA (pH 8.0) with blue dextran (50 mg/mL)</li></ul> <p><b>⚠ WARNING CHEMICAL HAZARD.</b> Formamide is harmful if absorbed through the skin and may cause irritation to the eyes, skin, and respiratory tract. It may cause damage to the central nervous system and the male and female reproductive systems, and is a possible birth defect hazard. Please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.</p> <p><b>⚠ CAUTION CHEMICAL HAZARD.</b> EDTA may cause eye, skin, and respiratory tract irritation. Please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.</p>
4	<p>Combine 4 <math>\mu\text{L}</math> of each 25-<math>\mu\text{L}</math> reaction/EDTA mixture with 4 <math>\mu\text{L}</math> of loading buffer.</p> <p>Store the remaining reaction/EDTA mixture at <math>-15</math> to <math>-25</math> <math>^{\circ}\text{C}</math>. This mixture can be concentrated by ethanol precipitation if the Express Load procedure does not yield enough signal.</p>
5	<p>Heat the samples at 98 <math>^{\circ}\text{C}</math> for 5 minutes with the lids open to denature and concentrate. Place on ice until ready to load.</p>
6	<p>Load 2.5 <math>\mu\text{L}</math> of each sample into a separate lane of the gel.</p>

## Electrophoresis on the 373 Instrument with BigDye Filter Wheel

**Requirements** General guidelines are provided below for running the ABI PRISM® 373 DNA Sequencers with the ABI PRISM® BigDye™ Filter Wheel<sup>2</sup> installed. For more detailed instructions, please refer the user's manual for your 373 instrument or to the user bulletin entitled *Using the ABI PRISM 373 BigDye Filter Wheel* (P/N 4304367).

### Gel

For 48-cm well-to-read (wtr), we recommended 5% Long Ranger™ gel.

### New Dye Set/Primer (Mobility) Files

Gel Formulation	Dye Set/Primer (Mobility) File
5% Long Ranger gel, 48-cm wtr <sup>a</sup>	DP373{BDv3-21M13}v1.mob
	DP373{BDv3-M13Rev}v1.mob

a. If you are running other wtr lengths, these are being tested. Please call Technical Support.

### Matrix Standards

**IMPORTANT** The instrument (matrix) file for the BigDye primers v3.0 cannot be used for the BigDye terminators (original), BigDye terminators v2.0, dRhodamine terminators, or BigDye primers (original).

Dye/Filter Set	Standards for Instrument (Matrix) File Generation
A (For use with the BigDye Filter Wheel)	ABI PRISM® BigDye™ Matrix Standards v3.0 (P/N 4390421)

**Note** For instructions on using the matrix standards (P/N 4390421) for the 373 instruments, contact Technical Support.

---

2. Includes the ABI PRISM 373 and ABI PRISM 373 with XL Upgrade instruments.

**Resuspending and Loading the Samples**

To resuspend and load the samples:

Step	Action																			
1	<p>Prepare a loading buffer by combining the following in a 5:1 ratio (5 parts deionized formamide to 1 part EDTA with blue dextran):</p> <ul style="list-style-type: none"> <li>• Deionized formamide</li> <li>• 25 mM EDTA (pH 8.0) with blue dextran (50 mg/mL)</li> </ul> <p><b>⚠ WARNING CHEMICAL HAZARD.</b> Formamide is harmful if absorbed through the skin and may cause irritation to the eyes, skin, and respiratory tract. It may cause damage to the central nervous system and the male and female reproductive systems, and is a possible birth defect hazard. Please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.</p> <p><b>⚠ CAUTION CHEMICAL HAZARD.</b> EDTA may cause eye, skin, and respiratory tract irritation. Please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.</p>																			
2	<p>Resuspend each sample pellet in loading buffer as follows:</p> <table border="1"> <thead> <tr> <th rowspan="2">Template</th> <th colspan="4">Volume (µL)</th> </tr> <tr> <th>18 or 24 well</th> <th>32 or 36 well</th> <th>48-well</th> <th>64-well</th> </tr> </thead> <tbody> <tr> <td>PCR product, plasmid, M13</td> <td>3–4</td> <td>3–4</td> <td>3–4</td> <td>2–4</td> </tr> <tr> <td>BAC, large DNA</td> <td>3</td> <td>3</td> <td>2</td> <td>2</td> </tr> </tbody> </table>	Template	Volume (µL)				18 or 24 well	32 or 36 well	48-well	64-well	PCR product, plasmid, M13	3–4	3–4	3–4	2–4	BAC, large DNA	3	3	2	2
Template	Volume (µL)																			
	18 or 24 well	32 or 36 well	48-well	64-well																
PCR product, plasmid, M13	3–4	3–4	3–4	2–4																
BAC, large DNA	3	3	2	2																
3	Vortex and spin the samples.																			
4	Heat the samples at 95 °C for 2 minutes to denature. Place on ice until ready to load.																			
5	<p>Load each sample into a separate lane of the gel as follows:</p> <table border="1"> <thead> <tr> <th rowspan="2">Template</th> <th colspan="4">Volume (µL)</th> </tr> <tr> <th>18 or 24 well</th> <th>32 or 36 well</th> <th>48-well</th> <th>64-well</th> </tr> </thead> <tbody> <tr> <td>PCR product, plasmid, M13</td> <td>1–3</td> <td>1–3</td> <td>1–2</td> <td>1–2</td> </tr> <tr> <td>BAC, large DNA</td> <td>3</td> <td>3</td> <td>2</td> <td>2</td> </tr> </tbody> </table>	Template	Volume (µL)				18 or 24 well	32 or 36 well	48-well	64-well	PCR product, plasmid, M13	1–3	1–3	1–2	1–2	BAC, large DNA	3	3	2	2
Template	Volume (µL)																			
	18 or 24 well	32 or 36 well	48-well	64-well																
PCR product, plasmid, M13	1–3	1–3	1–2	1–2																
BAC, large DNA	3	3	2	2																



# *Control DNA Sequences*



## **Appendix Summary**

---

**In This Appendix** The following topics are covered in this appendix:

<b>Topic</b>	<b>See Page</b>
-21 M13 Forward Primer	A-2
M13 Reverse Primer	A-3

---

## -21 M13 Forward Primer

**Sequence** The pGEM®-3Zf(+) sequence below is the sequence of the -21 M13 forward primer, followed by the ensuing 1000 bases.

---

TGTA AACGACGGCCAGT (-21 M13 primer)

GAATTGTAAT	ACGACTCACT	ATAGGGCGAA	TTCGAGCTCG	40
GTACCCGGGG	ATCCTCTAGA	GTCGACCTGC	AGGCATGCAA	80
GCTTGAGTAT	TCTATAGTGT	CACCTAAATA	GCTTGCGTA	120
ATCATGGTCA	TAGCTGTTTC	CTGTGTGAAA	TTGTTATCCG	160
CTCACAATTC	CACACAACAT	ACGAGCCGGA	AGCATAAAGT	200
GTA AAGCCTG	GGGTGCCTAA	TGAGTGAGCT	AACTCACATT	240
AATTGCGTTG	CGCTCACTGC	CCGCTTTCCA	GTCGGGAAAC	280
CTGTGCGGCC	AGCTGCATTA	ATGAATCGGC	CAACGCGCGG	320
GGAGAGGCGG	TTTGCGTATT	GGGCGCTCTT	CCGCTTCCTC	360
GCTCACTGAC	TCGCTGCGCT	CGGTCGTTCCG	GCTGCGGCGA	400
GCGGTATCAG	CTCACTCAA	GGCGGTAATA	CGGTTATCCA	440
CAGAATCAGG	GGATAACGCA	GGAAAGAACA	TGTGAGCAAA	480
AGGCCAGCAA	AAGGCCAGGA	ACCGTAAAAA	GGCCGCGTTG	520
CTGGCGTTTT	TCCATAGGCT	CCGCCCCCT	GACGAGCATC	560
ACAAAAATCG	ACGCTCAAGT	CAGAGGTGGC	GAAACCCGAC	600
AGGACTATAA	AGATACCAGG	CGTTTCCCCC	TGGAAGCTCC	640
CTCGTGCGCT	CTCCTGTTCC	GACCCTGCCG	CTTACCGGAT	680
ACCTGTCCGC	CTTTCTCCCT	TCGGGAAGCG	TGGCGCTTTC	720
TCATAGCTCA	CGCTGTAGGT	ATCTCAGTTC	GGTGTAGGTC	760
GTTGCTCCA	AGCTGGGCTG	TGTGCACGAA	CCCCCGTTC	800
AGCCCGACCG	CTGCGCCTTA	TCCGGA TACT	ATCGTCTTGA	840
GTCCAACCCG	GTAAGACACG	ACTTATCGCC	ACTGGCAGCA	880
GCCACTGGTA	ACAGGATTAG	CAGAGCGAGG	TATGTAGGCG	920
GTGCTACAGA	GTTCTTGAAG	TGGTGGCCTA	ACTACGGCTA	960
CACTAGAAGG	ACAGTATTTG	GTATCTGCGC	TCTGCTGAAG	1000

---



## M13 Reverse Primer

---

**Sequence** The pGEM®-3Zf(+) sequence below is the sequence of the M13 reverse primer, followed by the ensuing 1000 bases.

CAGGAAACAGCTATGACC (M13 reverse primer)

ATGATTACGC	CAAGCTATTT	AGGTGACACT	ATAGAATACT	40
CAAGCTTGCA	TGCCTGCAGG	TCGACTCTAG	AGGATCCCCG	80
GGTACCGAGC	TCGAATTCGC	CCTATAGTGA	GTCGTATTAC	120
AATTCACTGG	CCGTCGTTTT	ACAACGTCGT	GACTGGGAAA	160
ACCCTGGCGT	TACCCAACCTT	AATCGCCTTG	CAGCACATCC	200
CCCTTTCGCC	AGCTGGCGTA	ATAGCGAAGA	GGCCCGCACC	240
GATCGCCCTT	CCCAACAGTT	GCGCAGCCTG	AATGGCGAAT	280
GGACGCGCCC	TGTAGCGGCG	CATTAAGCGC	GGCGGGTGTG	320
GTGGTTACGC	GCAGCGTGAC	CGCTACACTT	GCCAGCGCCC	360
TAGCGCCCGC	TCCTTTCGCT	TTCTTCCCTT	CCTTTCTCGC	400
CACGTTTCGCC	GGCTTTCCCC	GTCAAGCTCT	AAATCGGGGG	440
CTCCCTTTAG	GGTTCGATT	TAGTGCTTTA	CGGCACCTCG	480
ACCCCAAAAA	ACTTGATTAG	GGTGATGGTT	CACGTAGTGG	520
GCCATCGCCC	TGATAGACGG	TTTTTCGCCC	TTTGACGTTG	560
GAGTCCACGT	TCTTTAATAG	TGGACTCTTG	TTCCAAACTG	600
GAACAACACT	CAACCCTATC	TCGGTCTATT	CTTTTGATTT	640
ATAAGGGATT	TTGCCGATTT	CGGCCTATTG	GTAAAAAAT	680
GAGCTGATTT	AACAAAAATT	TAACGCGAAT	TTTAACAAAA	720
TATTAACGCT	TACAATTTCC	TGATGCGGTA	TTTTCTCCTT	760
ACGCATCTGT	GCGGTATTTT	ACACCGCATA	TGGTGCACCT	800
TCAGTACAAT	CTGCTCTGAT	GCCGCATAGT	TAAGCCAGCC	840
CCGACACCCG	CCAACACCCG	CTGACGCGCC	CTGACGGGCT	880
TGTCTGCTCC	CGGCATCCGC	TTACAGACAA	GCTGTGACCG	920
TCTCCGGGAG	CTGCATGTGT	CAGAGGTTTT	CACCGTCATC	960
ACCGAAACGC	GCGAGACGAA	AGGGCCTCGT	GATACGCCTA	1000

---



# *Technical Support*

# B

## **Technical Support**

### **Contacting Technical Support**

---

You can contact Applied Biosystems for technical support by telephone or fax, by e-mail, or through the Internet. You can order Applied Biosystems user documents, MSDSs, certificates of analysis, and other related documents 24 hours a day. In addition, you can download documents in PDF format from the Applied Biosystems Web site (please see the section "To Obtain Documents on Demand" following the telephone information below).

---

### **To Contact Technical Support by E-Mail**

---

Contact technical support by e-mail for help in the following product areas:

<b>Product Area</b>	<b>E-mail address</b>
Genetic Analysis (DNA Sequencing)	galab@appliedbiosystems.com
Sequence Detection Systems and PCR	pcrlab@appliedbiosystems.com
Protein Sequencing, Peptide and DNA Synthesis	corelab@appliedbiosystems.com
Biochromatography, PerSeptive DNA, PNA and Peptide Synthesis systems, CytoFluor®, FMAT™, Voyager™, and Mariner™ Mass Spectrometers	tssupport@appliedbiosystems.com
Applied Biosystems/MDS Sciex	api3-support@sciex.com
Chemiluminescence (Tropix)	tropix@appliedbiosystems.com

---

**Hours for Telephone Technical Support**

In the United States and Canada, technical support is available at the following times:

Product	Hours
Chemiluminescence	8:30 a.m. to 5:30 p.m. Eastern Time
Framingham support	8:00 a.m. to 6:00 p.m. Eastern Time
All Other Products	5:30 a.m. to 5:00 p.m. Pacific Time

**To Contact Technical Support by Telephone or Fax**

**In North America**

To contact Applied Biosystems Technical Support, use the telephone or fax numbers given below. (To open a service call for other support needs, or in case of an emergency, dial **1-800-831-6844** and press **1**.)

Product or Product Area	Telephone Dial...	Fax Dial...
ABI PRISM® 3700 DNA Analyzer	<b>1-800-831-6844</b> , then press <b>8</b>	<b>1-650-638-5981</b>
DNA Synthesis	<b>1-800-831-6844</b> , then press <b>21</b>	<b>1-650-638-5981</b>
Fluorescent DNA Sequencing	<b>1-800-831-6844</b> , then press <b>22</b>	<b>1-650-638-5981</b>
Fluorescent Fragment Analysis (includes GeneScan® applications)	<b>1-800-831-6844</b> , then press <b>23</b>	<b>1-650-638-5981</b>
Integrated Thermal Cyclers (ABI PRISM®877 and Catalyst 800 instruments)	<b>1-800-831-6844</b> , then press <b>24</b>	<b>1-650-638-5981</b>
ABI PRISM® 3100 Genetic Analyzer	<b>1-800-831-6844</b> , then press <b>26</b>	<b>1-650-638-5981</b>
Bioinformatics (includes BioLIMS®, BioMerge®, and SQL GT™ applications)	<b>1-800-831-6844</b> , then press <b>25</b>	<b>1-505-982-7690</b>
Peptide Synthesis (433 and 43X Systems)	<b>1-800-831-6844</b> , then press <b>31</b>	<b>1-650-638-5981</b>
Protein Sequencing (Procise® Protein Sequencing Systems)	<b>1-800-831-6844</b> , then press <b>32</b>	<b>1-650-638-5981</b>

<b>Product or Product Area</b>	<b>Telephone Dial...</b>	<b>Fax Dial...</b>
PCR and Sequence Detection	<b>1-800-762-4001</b> , then press 1 for PCR, 2 for the 7700 or 5700, 6 for the 6700 or dial <b>1-800-831-6844</b> , then press 5	<b>1-240-453-4613</b>
Voyager™ MALDI-TOF Biospectrometry™ and Mariner™ ESI-TOF Mass Spectrometry Workstations	<b>1-800-899-5858</b> , then press 13	<b>1-508-383-7855</b>
Biochromatography (BioCAD® Workstations and POROS® Perfusion Chromatography Products)	<b>1-800-899-5858</b> , then press 14	<b>1-508-383-7855</b>
Expedite™ Nucleic acid Synthesis Systems	<b>1-800-899-5858</b> , then press 15	<b>1-508-383-7855</b>
Peptide Synthesis (Pioneer™ and 9050 Plus Peptide Synthesizers)	<b>1-800-899-5858</b> , then press 15	<b>1-508-383-7855</b>
PNA Custom and Synthesis	<b>1-800-899-5858</b> , then press 15	<b>1-508-383-7855</b>
FMAT™ 8100 HTS System and Cytofluor® 4000 Fluorescence Plate Reader	<b>1-800-899-5858</b> , then press 16	<b>1-508-383-7855</b>
Chemiluminescence (Tropix)	<b>1-800-542-2369</b> (U.S. only), or <b>1-781-271-0045</b>	<b>1-781-275-8581</b>
Applied Biosystems/MDS Sciex	<b>1-800-952-4716</b>	<b>1-650-638-6223</b>

### Outside North America

Region	Telephone Dial...	Fax Dial...
<b>Africa and the Middle East</b>		
Africa (English Speaking) and West Asia (Fairlands, South Africa)	27 11 478 0411	27 11 478 0349
South Africa (Johannesburg)	27 11 478 0411	27 11 478 0349
Middle Eastern Countries and North Africa (Monza, Italia)	39 (0)39 8389 481	39 (0)39 8389 493

<b>Eastern Asia, China, Oceania</b>		
Australia (Scoresby, Victoria)	61 3 9730 8600	61 3 9730 8799
China (Beijing)	86 10 64106608	86 10 64106617
Hong Kong	852 2756 6928	852 2756 6968
Korea (Seoul)	82 2 593 6470/6471	82 2 593 6472
Malaysia (Petaling Jaya)	60 3 758 8268	60 3 754 9043
Singapore	65 896 2168	65 896 2147
Taiwan (Taipei Hsien)	886 2 2358 2838	886 2 2358 2839
Thailand (Bangkok)	66 2 719 6405	66 2 319 9788
<b>Europe</b>		
Austria (Wien)	43 (0)1 867 35 75 0	43 (0)1 867 35 75 11
Belgium	32 (0)2 712 5555	32 (0)2 712 5516
Czech Republic and Slovakia (Praha)	420 2 61 222 164	420 2 61 222 168
Denmark (Naerum)	45 45 58 60 00	45 45 58 60 01
Finland (Espoo)	358 (0)9 251 24 250	358 (0)9 251 24 243
France (Paris)	33 (0)1 69 59 85 85	33 (0)1 69 59 85 00
Germany (Weiterstadt)	49 (0) 6150 101 0	49 (0) 6150 101 101
Hungary (Budapest)	36 (0)1 270 8398	36 (0)1 270 8288
Italy (Milano)	39 (0)39 83891	39 (0)39 838 9492
Norway (Oslo)	47 23 12 06 05	47 23 12 05 75

<b>Region</b>	<b>Telephone Dial...</b>	<b>Fax Dial...</b>
Poland, Lithuania, Latvia, and Estonia (Warszawa)	<b>48 (22) 866 40 10</b>	<b>48 (22) 866 40 20</b>
Portugal (Lisboa)	<b>351 (0)22 605 33 14</b>	<b>351 (0)22 605 33 15</b>
Russia (Moskva)	<b>7 095 935 8888</b>	<b>7 095 564 8787</b>
South East Europe (Zagreb, Croatia)	<b>385 1 34 91 927</b>	<b>385 1 34 91 840</b>
Spain (Tres Cantos)	<b>34 (0)91 806 1210</b>	<b>34 (0)91 806 1206</b>
Sweden (Stockholm)	<b>46 (0)8 619 4400</b>	<b>46 (0)8 619 4401</b>
Switzerland (Rotkreuz)	<b>41 (0)41 799 7777</b>	<b>41 (0)41 790 0676</b>
The Netherlands (Nieuwerkerk a/d IJssel)	<b>31 (0)180 331400</b>	<b>31 (0)180 331409</b>
United Kingdom (Warrington, Cheshire)	<b>44 (0)1925 825650</b>	<b>44 (0)1925 282502</b>
All other countries not listed (Warrington, UK)	<b>44 (0)1925 282481</b>	<b>44 (0)1925 282509</b>
<b>Japan</b>		
Japan (Hacchobori, Chuo-Ku, Tokyo)	<b>81 3 5566 6100</b>	<b>81 3 5566 6501</b>
<b>Latin America</b>		
Del.A. Obregon, Mexico	<b>305-670-4350</b>	<b>305-670-4349</b>

**To Reach  
Technical Support  
Through the  
Internet**

---

---

We strongly encourage you to visit our Web site for answers to frequently asked questions and for more information about our products. You can also order technical documents or an index of available documents and have them faxed or e-mailed to you through our site. The Applied Biosystems Web site address is

**<http://www.appliedbiosystems.com/techsupp>**

To submit technical questions from North America or Europe:

<b>Step</b>	<b>Action</b>
<b>1</b>	Access the Applied Biosystems Technical Support Web site.
<b>2</b>	Under the <b>Troubleshooting</b> heading, click <b>Support Request Forms</b> , then select the relevant support region for the product area of interest.
<b>3</b>	Enter the requested information and your question in the displayed form, then click <b>Ask Us RIGHT NOW</b> (blue button with yellow text).
<b>4</b>	Enter the required information in the next form (if you have not already done so), then click <b>Ask Us RIGHT NOW</b> .  You will receive an e-mail reply to your question from one of our technical experts within 24 to 48 hours.

---

---



**To Obtain Documents on Demand**

Free, 24-hour access to Applied Biosystems technical documents, including MSDSs, is available by fax or e-mail or by download from our Web site.

To order documents...	Then...
by index number	a. Access the Applied Biosystems Technical Support Web site at <b><a href="http://www.appliedbiosystems.com/techsupp">http://www.appliedbiosystems.com/techsupp</a></b> b. Click the <b>Index</b> link for the document type you want, then find the document you want and record the index number. c. Use the index number when requesting documents following the procedures below.
by phone for fax delivery	a. From the U.S. or Canada, call <b>1-800-487-6809</b> , or from outside the U.S. and Canada, call <b>1-858-712-0317</b> . b. Follow the voice instructions to order the documents you want. <b>Note</b> There is a limit of five documents per request.
through the Internet for fax or e-mail delivery	a. Access the Applied Biosystems Technical Support Web site at <b><a href="http://www.appliedbiosystems.com/techsupp">http://www.appliedbiosystems.com/techsupp</a></b> b. Under <b>Resource Libraries</b> , click the type of document you want. c. Enter or select the requested information in the displayed form, then click <b>Search</b> . d. In the displayed search results, select a check box for the method of delivery for each document that matches your criteria, then click <b>Deliver Selected Documents Now</b> (or click the PDF icon for the document to download it immediately). e. Fill in the information form (if you have not previously done so), then click <b>Deliver Selected Documents Now</b> to submit your order. <b>Note</b> There is a limit of five documents per request for fax deliver but no limit on the number of documents you can order for e-mail delivery.



## ABI PRISM® DNA Sequencing Kits and Related Products

To order ABI PRISM DNA sequencing kits, please contact Applied Biosystems at one of the regional sales offices listed on the back of this protocol. All reagents are quality-controlled in stable formulations. All the kits listed below include protocols. Protocols can also be ordered separately.

### dRhodamine Terminator Cycle Sequencing Kits with AmpliTaq® DNA Polymerase, FS

P/N	Kit	Reactions
403044	Ready Reaction	100
403045	Ready Reaction	1000
4303143	Ready Reaction	5000
403041	Protocol	–

### The ABI PRISM® BigDye™ Primer v3.0 Ready Reaction Cycle Sequencing Kit with AmpliTaq® DNA Polymerase, FS

P/N	Primer	Reactions
4390161	-21 M13	100
4390163	M13 reverse	100
4390157	-21 M13	5000
4390159	M13 reverse	5000

### The ABI PRISM® BigDye™ Terminator v3.0 Ready Reaction Cycle Sequencing Kit with AmpliTaq® DNA Polymerase, FS

P/N	Kit	Reactions
4390236	Ready Reaction	24
4390242	Ready Reaction	100
4390244	Ready Reaction	1000
4390246	Ready Reaction	5000
4390253	Ready Reaction	25000

### The ABI PRISM® dGTP BigDye™ Terminator v3.0 Ready Reaction Cycle Sequencing Kit with AmpliTaq® DNA Polymerase, FS

P/N	Kit	Reactions
4390229	Ready Reaction	100

### ABI PRISM® Lane Guide™ Lane Identification Kits for use on the 377 Sequencer

P/N	Kit	Reactions
4313682	Lane Guide	200
4313677	Lane Guide	1000
4313679	Lane Guide	5000
4313804	Protocol	–

### ABI PRISM Matrix Standards

P/N	Kit	Instrument
4390309	ABI PRISM® 3700 BigDye™ Terminator v3.0 Sequencing Standard	3700
4390303	ABI PRISM® BigDye™ Terminator v3.0 Sequencing Standard	3100/377/ 373/310
4390421	ABI PRISM® BigDye™ Matrix Standards v3.0	377/373/310



## Regional Sales Offices

### THE AMERICAS

United States  
Applied Biosystems  
850 Lincoln Centre Drive  
Foster City, CA 94404  
Tel: (650) 570-6667  
(800) 345-5224  
Fax: (650) 572-2743

Canada (Mississauga, Ontario)  
Tel: (905) 821-8183  
(800) 668-6913  
Fax: (905) 821-8246

Latin America (Del.A. Obregon,  
Mexico)  
Tel: (305) 670-4350  
Fax: (305) 670-4349

### EUROPE, AFRICA

Austria (Wien)  
Tel: 01 602 3101  
Fax: 01 602 5174

Benelux (Nieuwerkerk a/d IJssel,  
Netherlands)  
Tel: 31 (0)180-331400  
Fax: 31 (0)180-331409

Chechia Rep. (Praha)  
Tel: 2-61 22 21 64  
Fax: 2-61 22 21 68

Denmark (Allerød)  
Tel: 48 100 400  
Fax: 48 100 401

Finland (Espoo)  
Tel: 09 751 72 700  
Fax: 09 751 72 701

France (Paris)  
Tel: 33-1 69 59 85 85  
Fax: 33-1 69 59 85 00

Germany (Weiterstadt)  
Tel: (0) 6150/101-0  
Fax: (0) 6150/101-101

Hungary (Budapest)  
Tel: 36-1-258-8410  
Fax: 36-1-256-9802

Italy (Milano)  
Tel: (039) 23831  
Fax: (039) 2383492

Norway (Oslo)  
Tel: (0) 22 02 1500  
Fax: (0) 22 02 1501

Poland (Warszawa)  
Tel: (48 22) 866 4010  
Fax: (48 22) 866 4020

Portugal (Lisboa)  
Tel: (351-1) 386 0997  
Fax: (351-1) 386 1000

Russia (Moskva)  
Tel: 095 935 8888  
Fax: 095 564 8787

South Africa (Johannesburg)  
Tel: 27 11 478 0411  
Fax: 27 11 478 0349

Spain (Madrid)  
Tel: (91) 806 1200  
Fax: (91) 804 0414

Sweden (Stockholm)  
Tel: (0)8 619 4400  
Fax: (0)8 619 4401

Switzerland (Rotkreuz)  
Tel: 041-799 7708  
Fax: 041-790 0676

United Kingdom (Warrington,  
Cheshire)  
Tel: (01925) 825650  
Fax: (01925) 282502

### All Other European Countries, Middle East, West Asia, Africa Except South Africa

Langen, Germany  
Tel: 49 6103 708 301  
Fax: 49 6103 708 310

### EASTERN ASIA, CHINA, OCEANIA, PACIFIC RIM

Australia (Scoresby, Victoria)  
Tel: (03) 9212-8585  
Fax: (03) 9212-8502

China (Beijing)  
Tel: 86 10 6238 1156  
Fax: 86 10 6238 1162

Hong Kong  
Tel: 852 2756 6928  
Fax: 852 2756 6968

Japan (Chiba)  
Tel: (0473) 80-8500  
Fax: (0473) 80-8505

Korea (Seoul)  
Tel: 822 592 7238  
Fax: 822 532 4908

Malaysia (Kuala Lumpur)  
Tel: 60 3 758 1118  
Fax: 60 3 754 9043

Singapore  
Tel: 65 896 2118  
Fax: 65 896 2147

Taiwan (Taipei Hsish)  
Tel: 886 22 698 3505  
Fax: 886 22 698 3405

Thailand (Bangkok)  
Tel: 662 719 6406  
Fax: 662 319 9788