

PrepSEQ™ Rapid Spin Sample Preparation Kits

© Copyright 2009, Life Technologies Corporation. All rights reserved.

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

Information in this document is subject to change without notice. Life Technologies Corporation assumes no responsibility for any errors that may appear in this document. This document is believed to be complete and accurate at the time of publication. In no event shall Life Technologies Corporation be liable for incidental, special, multiple, or consequential damages in connection with or arising from the use of 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.

Limited Warranty for Consumables

Applied Biosystems warrants that the accompanying consumable product ("Consumable") will be free of defects in materials and workmanship upon delivery to buyer. All Consumable warranty claims must be made within one year from the date of delivery, but in any event no later than the earlier or lesser, as the case may be, of any shelf life date, expiration date, "use by" date, "guaranty date" or other end of recommended use date, or the number of uses, stated on the Consumable label or in Consumable literature accompanying the Consumable (the "Consumables Warranty Period"). Applied Biosystems will replace free of charge or refund the purchase price of any Consumable not meeting the above warranty if the warranty claim for the defective Consumable is received by Applied Biosystems within the Consumables Warranty Period. If the defects in material and workmanship diminished but did not prevent performance of the Consumable, any amount refunded may be apportioned according to Applied Biosystems' reasonable allocation of the value of the diminished performance. The decision of whether to replace the Consumable or refund the purchase price will be made by Applied Biosystems in its sole discretion. Notwithstanding the foregoing, custom Consumables made to specifications of the customer are sold "AS IS", without any warranty whatsoever, express or implied, except to the extent set forth in any separate express written limited warranty included in the documentation shipped with the Consumable or in Applied Biosystems' quotation. Except as set forth above, expiration dates, shelf life, "use by," guaranty or other end of recommended use dates are included for informational purposes only and shall not be deemed as a period of warranty. This limited warranty for Consumables is subject to the exclusions, conditions, exceptions, and limitations set forth below under the caption "EXCLUSIONS, CONDITIONS, EXCEPTIONS, AND LIMITATIONS."

EXCLUSIONS, CONDITIONS, EXCEPTIONS, AND LIMITATIONS

In addition to any exclusions, conditions, exceptions, or limitations set forth in any Applied Biosystems express written warranty statement, Applied Biosystems' warranties do not cover, and Applied Biosystems has no obligation replace or refund the purchase price of, any Consumable if a defect or failure arises or is caused by: Failure to store or use the product in accordance with Applied Biosystems' instructions; failure to use the product in accordance with good laboratory practices by technically qualified persons; other misuse or neglect, including but not limited to improper storage, maintenance, shipping, or handling, improper or abnormal use, including operation outside of the environmental or use specifications specified by Applied Biosystems; adulteration; unauthorized change or modification; use with chemicals, reagents or any other product not supplied or authorized in writing by Applied Biosystems for use with the product; and accidents of any kind or nature.

Buyer is responsible for any validation of assays and compliance with any regulatory requirements that pertain to their procedures and uses of the Products. Improper or unapproved or noncompliant uses of the Products render any Applied Biosystems obligations under this warranty null and void.

Any claims related to breach of warranty not made by the customer within the applicable warranty period will be forfeited. Warranties are limited to the buyer of the product from Applied Biosystems and are not transferable.

THE FOREGOING WARRANTIES ARE APPLIED BIOSYSTEMS' SOLE AND EXCLUSIVE WARRANTIES WITH RESPECT TO INSTRUMENTS, COMPONENTS, SOFTWARE, SERVICE PARTS AND CONSUMABLES, AND IS IN LIEU OF ALL OTHER WARRANTIES, EXPRESSED OR IMPLIED, ALL OF WHICH OTHER WARRANTIES ARE EXPRESSLY DISCLAIMED, INCLUDING WITHOUT LIMITATION THOSE OF MERCHANTABILITY, SATISFACTORY QUALITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT, WHETHER ARISING FROM A STATUTE OR OTHERWISE IN LAW OR FROM A COURSE OF DEALING OR USAGE OF TRADE. REPAIR, REPLACEMENT OR REFUND OF THE PURCHASE PRICE AS SET FORTH ABOVE IS APPLIED BIOSYSTEMS' SOLE OBLIGATION AND CUSTOMER'S SOLE AND EXCLUSIVE REMEDY WITH RESPECT TO APPLIED BIOSYSTEMS' BREACH OF WARRANTY.

IN NO EVENT SHALL APPLIED BIOSYSTEMS BE LIABLE, WHETHER IN CONTRACT, TORT, STRICT LIABILITY, NEGLIGENCE, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR:

SPECIAL LOSSES OR DAMAGES;

INCIDENTAL LOSSES OR DAMAGES;

INDIRECT LOSSES OR DAMAGES;

EXEMPLARY, PUNITIVE, OR MULTIPLE LOSSES OR DAMAGES;

CONSEQUENTIAL LOSSES OR DAMAGES, OR

INCREASED COSTS OR EXPENSES

SUSTAINED BY CUSTOMER OR ANY OTHER PERSON OR ENTITY ARISING OUT OF OR CAUSED BY THE PURCHASE, POSSESSION OR USE OF APPLIED BIOSYSTEMS' PRODUCTS OR PRODUCTS SUPPLIED BY APPLIED BIOSYSTEMS, OR THE PERFORMANCE BY APPLIED BIOSYSTEMS OF ANY SERVICES RELATING TO ANY SUCH PRODUCTS, WHETHER OR NOT FORESEEABLE AND WHETHER

OR NOT APPLIED BIOSYSTEMS IS ADVISED OF THE POSSIBILITY OF SUCH DAMAGES, INCLUDING WITHOUT LIMITATION DAMAGES ARISING FROM OR RELATED TO LOSS OF USE, LOSS OF DATA, DOWNTIME, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES, OR FOR LOSS OF REVENUE, PROFITS, INCOME, GOODWILL, OR BUSINESS OR OTHER FINANCIAL LOSS OR PERSONAL INJURY OR PROPERTY DAMAGE, INFRINGEMENT, OR REGARDING RESULTS OBTAINED THROUGH THE USE OF ANY PRODUCT (INCLUDING WITHOUT LIMITATION, ANY CLAIM OF INACCURATE, INVALID, OR INCOMPLETE RESULTS).

Some countries or jurisdictions limit the scope of or preclude limitations or exclusion of warranties, of liability, such as liability for gross negligence or willful misconduct, or of remedies or damages, as or to the extent set forth above. In such countries and jurisdictions, the limitation or exclusion of warranties, liability, remedies or damages set forth above shall apply to the fullest extent permitted by law, and shall not apply to the extent prohibited by law. Without limiting the foregoing statement, nothing in this document is intended to or will exclude or limit Applied Biosystems' liability for death or personal injury caused by its negligence or fraud, or any other liability to the extent that the same may not be excluded or limited by applicable law.

NOTICE TO PURCHASER: License Disclaimer

This product conveys no patent rights, expressly or by implication, under any patent or patent application owned by or licensable by Applied Biosystems that covers any thermal cycling instrument, apparatus or system, any composition, reagent, or kit, or any process. Specifically, but without limitation, no right, immunity, authorization, or license is granted, expressly or by implication, for the processes of PCR, real-time PCR, reverse-transcription PCR, or the 5' nuclease assay.

Further information regarding 5' nuclease licensing program and commercial service licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

TRADEMARKS:

The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

	Preface	v
	Safety information	v
	How to use this guide	vi
	How to obtain support	vi
Protocol	PrepSEQ™ Rapid Spin Sample Preparation Kits	1
	Product overview	1
	Required materials	1
	Food sample preparation	4
	Troubleshooting	14
Appendix A	Safety	15
	Chemical safety	16
	Chemical waste safety	17
	Biological hazard safety	19
	Chemical alerts	19
	Documentation	21
	Related documentation	21

Safety information

Note: For general safety information, see this Preface and [Appendix A, “Safety” on page 15](#). When a hazard symbol and hazard type appear by a chemical name or instrument hazard, see the “Safety” Appendix for the complete alert on the chemical or instrument.

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:

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.

MSDSs

The MSDSs for any chemicals supplied by Applied Biosystems or Ambion are available to you free 24 hours a day. For instructions on obtaining MSDSs, see [Appendix A](#).

IMPORTANT! For the MSDSs of chemicals not distributed by Applied Biosystems or Ambion contact the chemical manufacturer.

How to use this guide

Text conventions

This guide uses the following conventions:

- **Bold** text indicates user action. For example:
Type **0**, then press **Enter** for each of the remaining fields.
- *Italic* text indicates new or important words and is also used for emphasis.
For example:
Before analyzing, *always* prepare fresh matrix.
- A right arrow symbol (▶) separates successive commands you select from a drop-down or shortcut menu. For example:
Select **File ▶ Open ▶ Spot Set**.
Right-click the sample row, then select **View Filter ▶ View All Runs**.

User attention words

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

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

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

How to obtain support

For the latest services and support information for all locations, go to www.appliedbiosystems.com.

At the Applied Biosystems web site, you can:

- Access 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

PrepSEQ™ Rapid Spin Sample Preparation Kits

Product overview

The PrepSEQ™ Rapid Spin Sample Preparation Kits provide a simple way to prepare DNA from broth cultures.

Required materials

Kit contents

The PrepSEQ™ Rapid Spin Sample Preparation Kits contain reagents for 100 sample preparations. Kit components are shown in the table below. For information on the kit contents, refer to the “Materials supplied” section in the packaging insert for your kit.

Kit	Item	Quantity or volume
Kits without Proteinase K		
PrepSEQ™ Rapid Spin Sample Preparation Kit (PN 4407760)	Spin columns	100
	Microcentrifuge tubes, 1.5 mL	100
	Lysis Buffer, 1 bottle	5 mL
PrepSEQ™ Rapid Spin Sample Preparation Kit – Extra Clean (PN 4413269)	Spin columns	100
	Microcentrifuge tubes, 1.5 mL	2 × 100
	Lysis Buffer, 1 bottle	5 mL
Kits with Proteinase K for difficult-to-lyse bacteria[‡]		
PrepSEQ™ Rapid Spin Sample Preparation Kit with Proteinase K (PN 4426714)	Spin columns	100
	Microcentrifuge tubes, 1.5 mL	100
	Lysis Buffer, 1 bottle	5 mL
	Proteinase K [‡] (20 mg/mL), 1 tube	1.25 mL
PrepSEQ™ Rapid Spin Sample Preparation Kit – Extra Clean with Proteinase K (PN 4426715)	Spin columns	100
	Microcentrifuge tubes, 1.5 mL	2 × 100
	Lysis Buffer, 1 bottle	5 mL
	Proteinase K [‡] (20 mg/mL), 1 tube	1.25 mL

[‡] for use with the protocol for difficult-to-lyse bacteria, such as *Listeria* spp.

Note: Parts may ship separately depending on the configuration ordered and storage conditions.

Storage

- Store Proteinase K at –20 °C.
- Store Lysis Buffer at 4 °C.
- Store spin columns and tubes at room temperature (22 to 28 °C).

For information on storage of individual kit components, refer to the “Storage” section in the packaging insert for your kit.

Materials not included in the kit

The following table includes materials and equipment for using (but not included in) the PrepSEQ™ Rapid Spin Sample Preparation Kits. Unless otherwise indicated, many of the listed items are available from major laboratory suppliers (MLS).

Equipment, consumables, and reagents	
Item	Source
Equipment	
Block heater, 95 °C	MLS
Block heater, 56 °C [‡]	MLS
Rack for 1.5-mL tubes	MLS
Benchtop microcentrifuge (Eppendorf 5415 D or equivalent)	MLS
Stomacher 400 Laboratory Blender	Seward #0400/001/AJ
Vortexer	MLS
Consumables	
Disposable gloves	MLS
Micropipette tips, aerosol-resistant	MLS
Pipettors: <ul style="list-style-type: none"> • Positive-displacement • Air-displacement 	MLS
Whirl-Pak Filter Bags, 6" × 9", 24 oz, 250/pkg (Stomacher bags with mesh)	VWR #11216-520
Whirl-Pak Filter Bags, 6" × 9", 24 oz (Stomacher bags without mesh)	VWR #11216-280
Reagents	
Bacterial growth media [§]	MLS

Equipment, consumables, and reagents (continued)

Item	Source
Nuclease-free Water	Applied Biosystems PN AM9938

‡ for use with protocol for difficult-to-lyse bacteria, such as *Listeria* spp.

§ microorganism appropriate

Food sample preparation

Overview

The PrepSEQ™ Rapid Spin Sample Preparation Kits are designed to work with most food types. The kit procedure involves:

- Sample enrichment
- Sample preparation

For sample preparation from enriched food samples, Applied Biosystems recommends a 750-µL sample volume.

For the detection of most food-borne pathogens, use the “[General sample preparation protocol](#)” on page 8 with either the PrepSEQ™ Rapid Spin Sample Preparation Kit (PN 4407760) or PrepSEQ™ Rapid Spin Sample Preparation Kit – Extra Clean (PN 4413269).

For the detection of difficult-to-lyse bacteria (for example, *Listeria* spp.), use the “[Sample preparation protocol for difficult-to-lyse bacteria](#)” on page 11 with either the PrepSEQ™ Rapid Spin Sample Preparation Kit with Proteinase K (PN 4426714) or PrepSEQ™ Rapid Spin Sample Preparation Kit – Extra Clean with Proteinase K (PN 4426715).

IMPORTANT! Proteinase K is required for efficient lysis of *Listeria monocytogenes*.

For some foods with a high lipid content, such as infant formula, soft cheese, whole milk, smoked salmon (lox), and chicken wing samples, use the PrepSEQ™ Rapid Spin Extra Clean protocol with either the PrepSEQ™ Rapid Spin Sample Preparation Kit – Extra Clean (PN 4413269) or PrepSEQ™ Rapid Spin Sample Preparation Kit – Extra Clean with Proteinase K (PN 4426715).

For information on preparation of:

- *Listeria monocytogenes* DNA from food samples, refer to the *PrepSEQ™ Rapid Spin Sample Preparation Kit Protocol: Listeria monocytogenes* (PN 4412851)
- *Salmonella* spp. DNA from food samples, refer to the *PrepSEQ™ Rapid Spin Sample Preparation Kit Protocol: Salmonella* spp. (PN 4412848)

Before you begin

Before starting your sample extraction:

- If you are using the general sample preparation protocol:
 - Set the block heater temperature to 95 °C.
 - Label 1.5-mL microcentrifuge tubes.
- If you are using the sample preparation protocol for difficult-to-lyse bacteria:
 - Set one block heater temperature to 95 °C, and the other block heater temperature to 56 °C.
 - Label 1.5-mL microcentrifuge tubes.

- Prepare Proteinase K-Lysis Buffer mix: premix 5 µL of Proteinase K (20 mg/mL) with 50 µL of Lysis Buffer for each sample (use a clean appropriately-sized container for mixing). Multiply volumes by the number of samples plus 10% for overage. Mix well to disperse Proteinase K in Lysis Buffer. Store on ice until ready to use.

Kit workflow using the general sample preparation protocol

The figure below shows a sample processing workflow based on the general sample preparation protocol. For details, see [page 8](#).

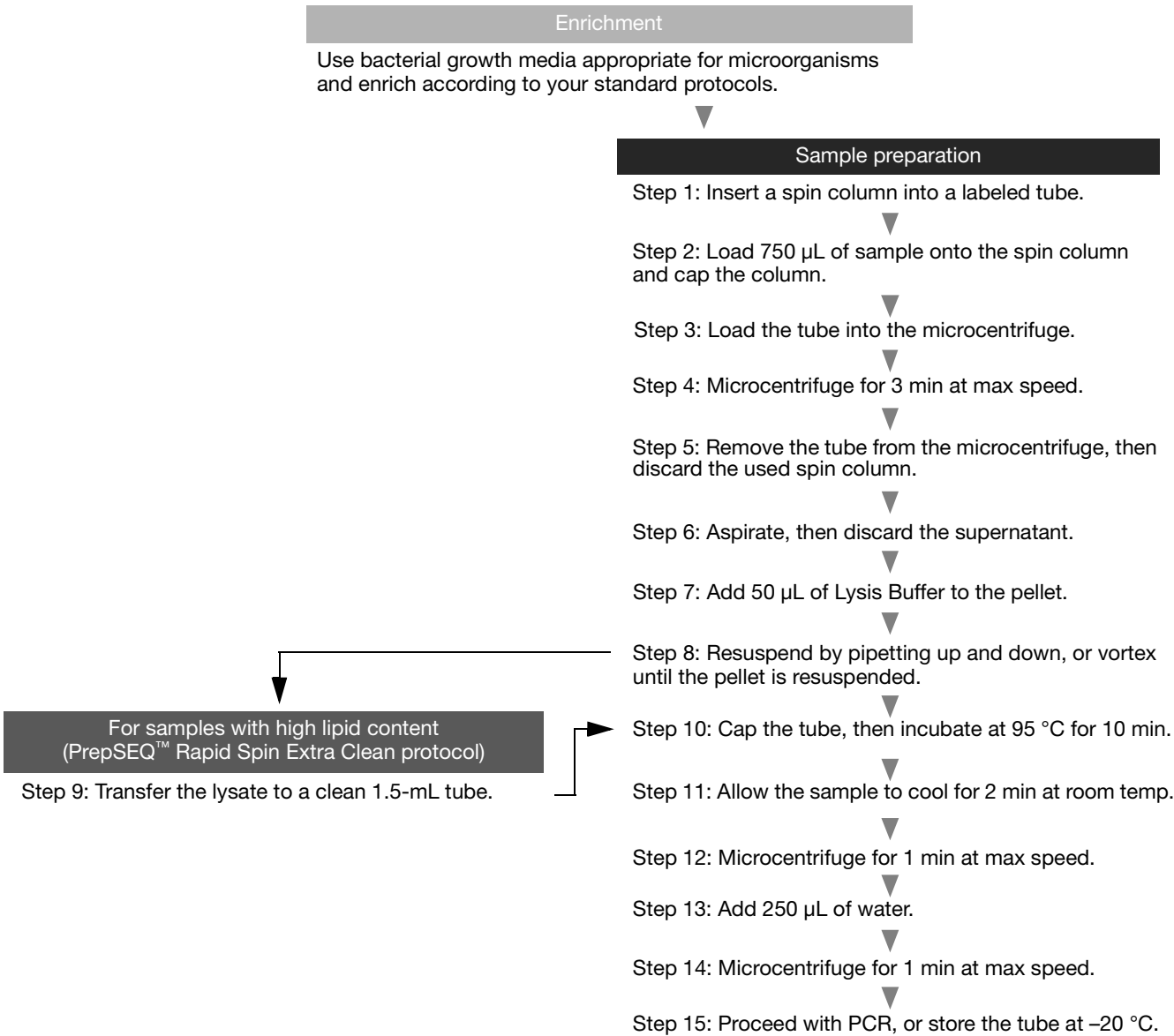


Figure 1 Food sample preparation using the PrepSEQ™ Rapid Spin Sample Preparation Kit general protocol

Kit workflow using the sample preparation protocol for difficult-to-lyse bacteria

The figure below shows a sample processing workflow based on the sample preparation protocol for difficult-to-lyse bacteria, such as *Listeria* spp. For details, see [page 11](#).

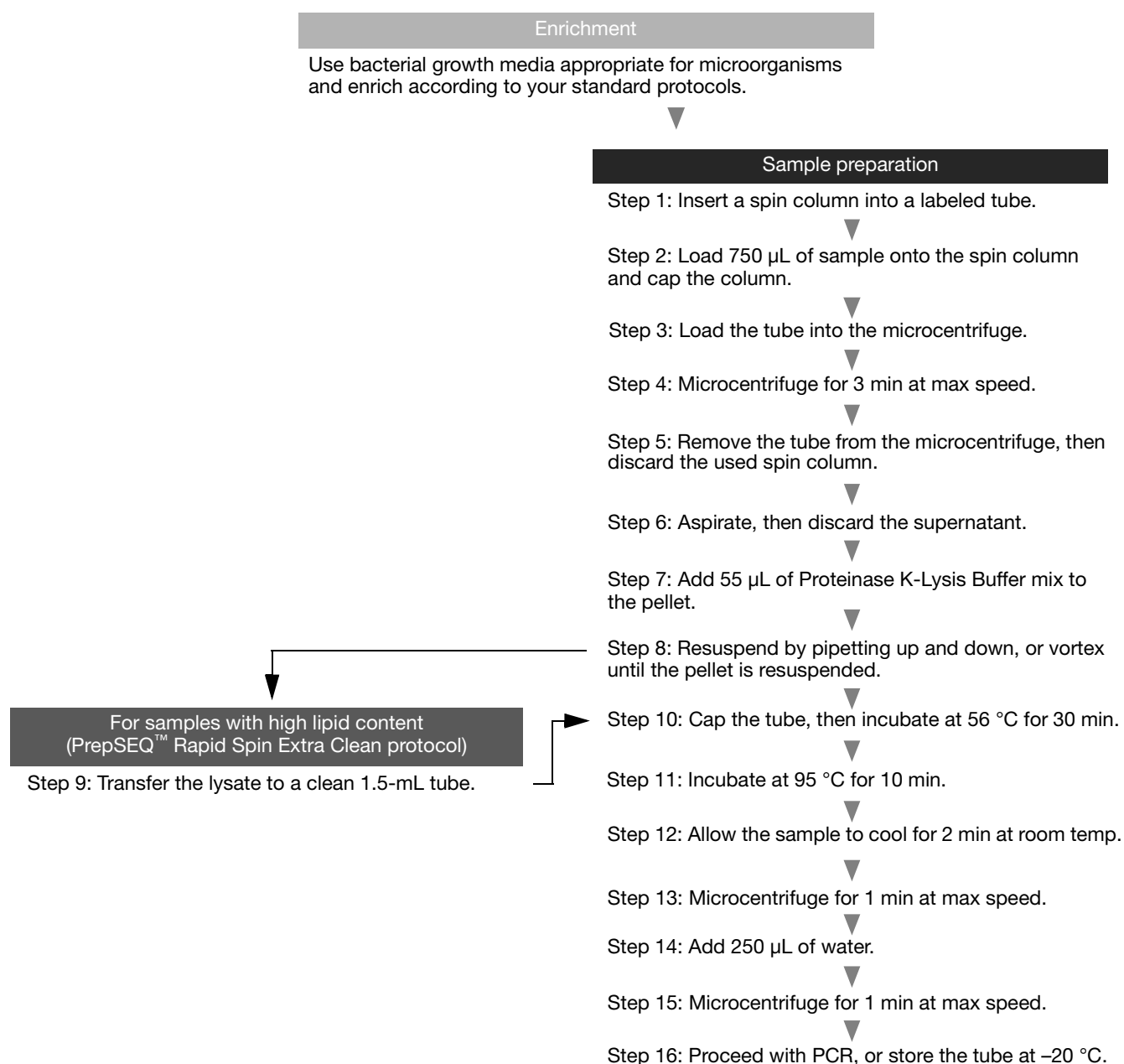


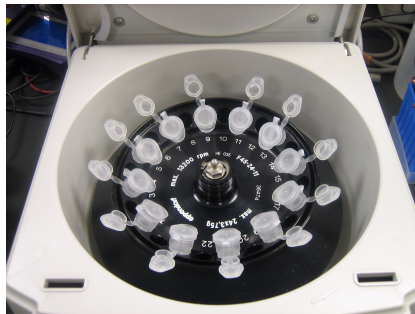
Figure 2 Food sample preparation using the PrepSEQ™ Rapid Spin Sample Preparation Kit protocol for difficult-to-lyse bacteria

Generic enrichment

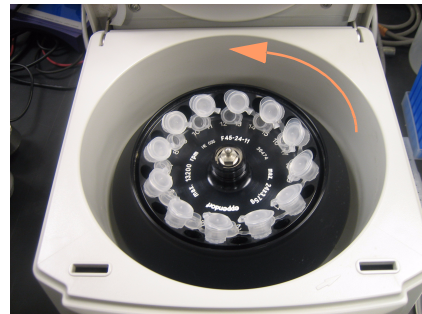
Use bacterial growth media appropriate for microorganisms and enrich according to your standard protocols.

General sample preparation protocol

1. Insert a spin column into a labeled tube.
2. Load 750 µL of your enriched sample onto the spin column and cap the column.
3. Load the tube with its spin column into the microcentrifuge. Place the tube cap hinge toward the inside of the rotor (see figure at right below), otherwise the cap hinge interferes with installation of the rotor lid.



Incorrect position of the tube cap



Correct position of the tube cap

4. Microcentrifuge the tube for 3 minutes at maximum speed.
5. Remove the tube from the microcentrifuge and discard the used spin column.
6. Aspirate, then discard the supernatant.

IMPORTANT! Remove supernatant as completely as possible, including any excess liquid on the sides of the tube. Remove droplets by circling the inside of the tube with the pipettor and pushing into supernatant for removal by aspiration.

IMPORTANT! For samples that contain a fat layer following centrifugation, indicated as a distinct top layer, remove the fat layer as follows:

- For liquid fat layer (for example, as found in soft cheese samples), use a pipettor to remove fat from the top surface by aspirating in a circular motion. Continue to collect supernatant from the top surface until all the supernatant is removed (discard into a waste container).
 - For solid fat layer (for example, as found in infant formula samples), use a pipettor to gently dislodge the fat layer and pour off the supernatant and fat layer using a quick motion (discard into a waste container). Remove the remaining supernatant using a pipettor.
-

7. Add 50 µL of Lysis Buffer to the pellet.

8. Resuspend by pipetting up and down, or vortex until the pellet is well dispersed in the Lysis Buffer.
9. Rapid Spin Extra Clean protocol *only* – If the food sample has high lipid content, transfer the Lysis Buffer mixture into a clean 1.5-mL tube (avoid transferring the fat when transferring the mixture). The pellet must be well dispersed in the Lysis Buffer prior to transfer. For all other samples, proceed directly to [step 10](#).

IMPORTANT! During the transfer there will be residual fat on the sides of the original tube. Avoid contact with the fat and transfer only the Lysis Buffer containing the resuspended pellet into a clean tube.

IMPORTANT! Applied Biosystems recommends this extra clean step for food samples with high lipid content, such as infant formula, soft cheese, whole milk, smoked salmon (lox), and chicken wing samples; for use with the PrepSEQ™ Rapid Spin Sample Preparation Kit – Extra Clean with Proteinase K (PN 4426715).

10. Cap the tube, then incubate at 95 °C for 10 minutes.
11. Allow the sample to cool for 2 minutes at room temperature.
12. Microcentrifuge the tube for 1 minute at maximum speed in order to bring down condensation following the 95 °C heating step.
13. Add 250 µL of water. Mix well.

IMPORTANT! Use PCR-clean water (PN AM9938). Autoclaved water should not be considered PCR-clean.

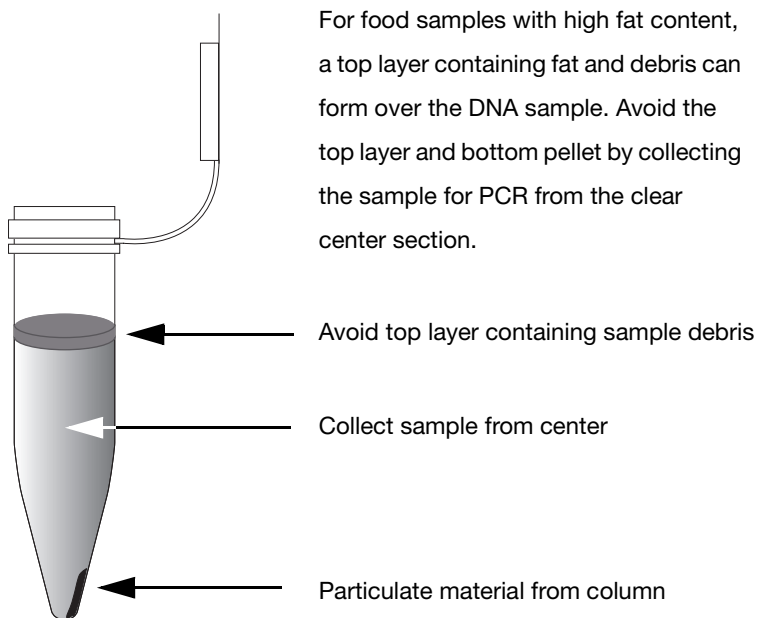
14. Microcentrifuge the tube for 1 minute at maximum speed to bring down any particulate material derived from the spin column, which can interfere with amplification. The microbial DNA is in the aqueous phase.

15. Proceed with PCR, or store the tube at -20°C .

IMPORTANT! Use 30 μL of supernatant for PCR in the lyophilized assay.

IMPORTANT! Food samples with high fat content can form a top layer containing fat and debris over the DNA sample. When collecting your sample for PCR avoid the top layer and bottom pellet by collecting your sample from the clear center section (see figure below).

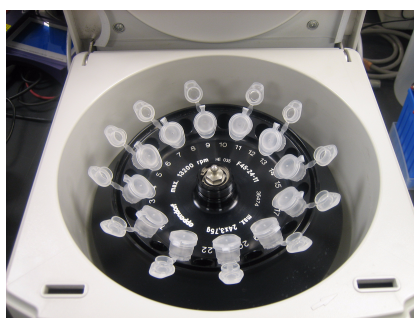
IMPORTANT! If PCR inhibition occurs, as indicated by no amplification for either target or IPC, then dilute sample with water. It is recommended to add 5 μL of sample and 25 μL of water to lyophilized assay to overcome inhibition. See [“Troubleshooting” on page 14](#) for further discussion.



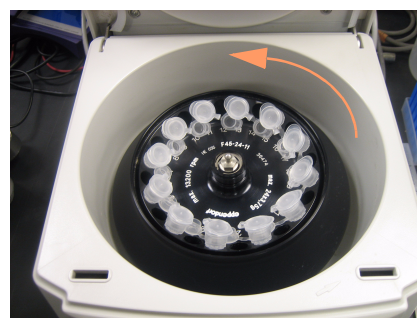
Sample preparation protocol for difficult-to-lyse bacteria

The following sample preparation protocol is for difficult-to-lyse bacteria, such as *Listeria* spp.

1. Insert a spin column into a labeled tube.
2. Load 750 µL of your enriched sample onto the spin column and cap the column.
3. Load the tube with its spin column into the microcentrifuge. Place the tube cap hinge toward the inside of the rotor (see figure at right below), otherwise the cap hinge interferes with installation of the rotor lid.



Incorrect position of the tube cap



Correct position of the tube cap

4. Microcentrifuge the tube for 3 minutes at maximum speed.
5. Remove the tube from the microcentrifuge and discard the used spin column.
6. Aspirate, then discard the supernatant.

IMPORTANT! Remove supernatant as completely as possible, including any excess liquid on the sides of the tube. Remove droplets by circling the inside of the tube with the pipettor and pushing into supernatant for removal by aspiration.

IMPORTANT! For samples that contain a fat layer following centrifugation, indicated as a distinct top layer, remove the fat layer as follows:

- For liquid fat layer (for example, as found in soft cheese samples), use a pipettor to remove fat from the top surface by aspirating in a circular motion. Continue to collect supernatant from the top surface until all the supernatant is removed (discard into a waste container).
 - For solid fat layer (for example, as found in infant formula samples), use a pipettor to gently dislodge the fat layer and pour off the supernatant and fat layer using a quick motion (discard into a waste container). Use a new pipette tip to remove the remaining supernatant.
-

7. Add 55 µL of Proteinase K-Lysis Buffer mix to the pellet.

IMPORTANT! To prepare the Proteinase K-Lysis Buffer mix: premix 5 µL of Proteinase K (20 mg/mL) with 50 µL of Lysis Buffer for each sample (use a clean appropriately-sized container for mixing). Multiply volumes by the number of samples plus 10% for overage. Mix well to disperse Proteinase K in Lysis Buffer.

IMPORTANT! Store Proteinase K-Lysis Buffer mix on ice until ready to use.

IMPORTANT! Proteinase K is required for efficient lysis of *Listeria* spp.

8. Resuspend by pipetting up and down, or vortex until the pellet is well dispersed in the Proteinase K-Lysis Buffer mix.
9. Rapid Spin Extra Clean with Proteinase K protocol *only* – If the food sample has high lipid content, transfer the mixture into a clean 1.5-mL tube (avoid transferring the fat when transferring the mixture). The pellet must be well dispersed in the Lysis Buffer prior to transfer.
For all other samples, proceed directly to [step 10](#) below.

IMPORTANT! During the transfer there will be residual fat on the sides of the original tube. Avoid contact with the fat and transfer only the Lysis Buffer containing the resuspended pellet into a clean tube.

IMPORTANT! Applied Biosystems recommends this extra clean step for food samples with high lipid content, such as infant formula, soft cheese, whole milk, smoked salmon (lox), and chicken wing samples; for use with the PrepSEQ™ Rapid Spin Sample Preparation Kit – Extra Clean with Proteinase K (PN 4426715).

10. Cap the tube, then incubate at 56 °C for 30 minutes.
11. Incubate at 95 °C for 10 minutes.
12. Allow the sample to cool for 2 minutes at room temperature.
13. Microcentrifuge the tube for 1 minute at maximum speed in order to bring down condensation following the 95 °C heating step.
14. Add 250 µL of water. Mix well.

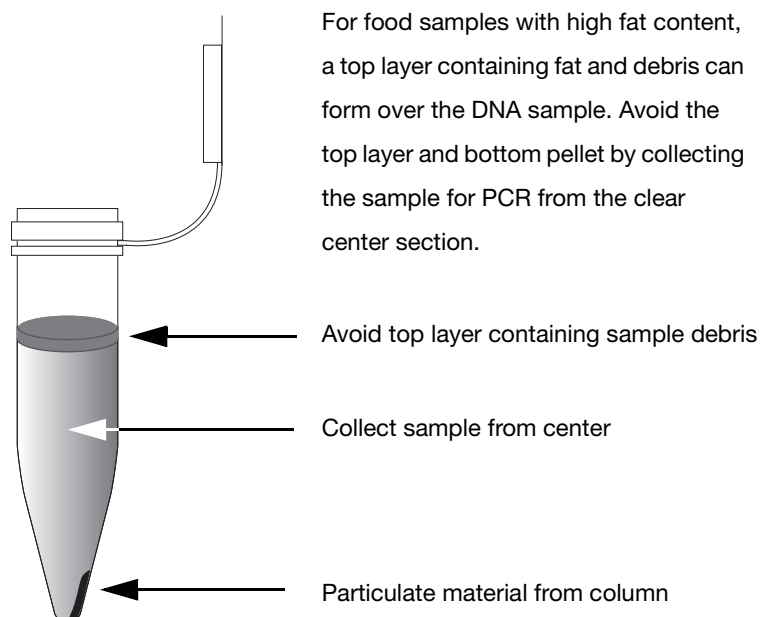
IMPORTANT! Use PCR-clean water (PN AM9938). Autoclaved water should not be considered PCR-clean.

15. Microcentrifuge the tube for 1 minute at maximum speed to bring down any particulate material derived from the spin column, which can interfere with amplification. The microbial DNA is in the aqueous phase.
16. Proceed with PCR, or store the tube at -20°C .

IMPORTANT! Use 30 μL of supernatant for PCR in the lyophilized assay.

IMPORTANT! Food samples with high fat content can form a top layer containing fat and debris over the DNA sample. When collecting your sample for PCR avoid the top layer and bottom pellet by collecting your sample from the clear center section (see figure below).

IMPORTANT! If PCR inhibition occurs, as indicated by no amplification for either target or IPC, then dilute sample with water. It is recommended to add 5 μL of sample and 25 μL of water to lyophilized assay to overcome inhibition. See [“Troubleshooting” on page 14](#) for further discussion.



Troubleshooting

Observation	Possible cause	Action
PCR is inhibited, indicated by non-detection of IPC reaction	Removal of the supernatant before adding Lysis Buffer was not sufficient.	Add 5 µL or 10 µL of sample to the PCR reaction and bring the final volume to 30 µL with water.
	Filtrate from the spin column is in the sample.	Centrifuge the sample to separate the filter particulates before adding sample to the PCR reaction.
	The sample contained excess fat that was not removed during aspiration of the supernatant.	Apply PrepSEQ™ Rapid Spin Extra Clean protocol.
	Matrix is associated with PCR inhibitory components.	Pre-wash the bacterial pellet before column clarification: 1 – Transfer 750 µL of sample to a clean microcentrifuge tube. 2 – Centrifuge at 16000 × g for 3 min. 3 – Discard supernatant. 4 – Resuspend pellet in 650 µL of sterile distilled water. 5 – Load column.
The bacterial pellet separates from the tube making pellet hard to avoid during aspiration	Sample was left unattended before the supernatant was aspirated.	Remove supernatant immediately following centrifugation.
Positive signal in negative control	Target organism is causing contamination.	PCR is a very sensitive test capable of amplification from a single copy of DNA. PCR must be set up in a PCR-clean environment that is separate from PCR amplification and analysis.

Safety

This appendix covers:

■ Chemical safety.....	16
■ Chemical waste safety.....	17
■ Biological hazard safety	19
■ Chemical alerts	19



Chemical safety

Chemical hazard warning



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



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



WARNING! CHEMICAL HAZARD. Four-liter reagent and waste bottles can crack and leak. Each 4-liter 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.



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

Chemical safety guidelines

To minimize the hazards of chemicals:

- Read and understand the Material Safety Data Sheets (MSDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. (See [“About MSDSs” on page 17.](#))
- 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 in the MSDS.
- Comply with all local, state/provincial, or national laws and regulations related to chemical storage, handling, and disposal.

About MSDSs

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

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

Obtaining MSDSs

The MSDS for any chemical supplied by Applied Biosystems is available to you free 24 hours a day. To obtain MSDSs:

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

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

Chemical waste safety

Chemical waste hazards



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



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



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



Chemical waste safety guidelines

To minimize the hazards of chemical waste:

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

Waste disposal

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

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

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

Biological hazard safety

General biohazard



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

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

Additional information about biohazard guidelines is available at: www.cdc.gov

Chemical alerts

General alerts for all chemicals

Avoid contact with skin, eyes, and/or clothing. Read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.



Related documentation

For additional documentation, see “How to obtain support” on page vi.

For information on new assays and updated product documentation, go to <http://info.appliedbiosystems.com/pathogenkits>

Document title	PN
<i>PrepSEQ™ Rapid Spin Sample Preparation Kit Quick Reference Card</i>	4412846
<i>PrepSEQ™ Rapid Spin Sample Preparation Kit Protocol: Salmonella spp.</i>	4412848
<i>PrepSEQ™ Rapid Spin Sample Preparation Kit Quick Reference Card: Salmonella spp.</i>	4412849
<i>PrepSEQ™ Rapid Spin Sample Preparation Kit Protocol: Listeria monocytogenes</i>	4412851
<i>PrepSEQ™ Rapid Spin Sample Preparation Kit Quick Reference Card: Listeria monocytogenes</i>	4412852
<i>PrepSEQ™ Nucleic Acid Extraction Kit Quick Reference Card</i>	4406303
<i>PrepSEQ™ Nucleic Acid Extraction Kit Protocol: Salmonella spp.</i>	4405968
<i>PrepSEQ™ Nucleic Acid Extraction Kit Quick Reference Card: Salmonella spp.</i>	4405967
<i>PrepSEQ™ Nucleic Acid Extraction Kit Protocol: Listeria monocytogenes</i>	4405966
<i>PrepSEQ™ Nucleic Acid Extraction Kit Quick Reference Card: Listeria monocytogenes</i>	4405965
<i>PrepSEQ™ Mycoplasma Nucleic Acid Extraction Kit Protocol</i>	4401253
<i>PrepSEQ™ Mycoplasma Nucleic Acid Extraction Kit Quick Reference Card</i>	4406304
<i>MicroSEQ® Listeria monocytogenes Detection Kit Protocol</i>	4405962
<i>MicroSEQ® Listeria monocytogenes Detection Kit Quick Reference Card</i>	4405961
<i>MicroSEQ® Salmonella spp. Detection Kit Protocol</i>	4405964
<i>MicroSEQ® Salmonella spp. Detection Kit Quick Reference Card</i>	4405963
<i>MycoSEQ™ Mycoplasma Detection Kits Protocol: MycoSEQ™ Mycoplasma Real-Time PCR Detection Kit, MycoSEQ™ Myco Scan Mycoplasma Detection Kit</i>	4393111
<i>MycoSEQ™ Mycoplasma Detection Kits Quick Reference Card: MycoSEQ™ Mycoplasma Real-Time PCR Detection Kit, MycoSEQ™ Myco Scan Mycoplasma Detection Kit</i>	4393471

Portable document format (PDF) versions of this guide and the documents listed above are available at www.appliedbiosystems.com

Note: To open the documentation available from the Applied Biosystems web site, use the Adobe® Acrobat® Reader® software available at www.adobe.com

Headquarters

850 Lincoln Centre Drive | Foster City, CA 94404 USA
Phone 650.638.5800 | Toll Free 800.345.5224
www.appliedbiosystems.com

International Sales

For our office locations please call the division
headquarters or refer to our Web site at
www.appliedbiosystems.com/about/offices.cfm