

PrepSEQ™ Rapid Spin Sample Preparation Kit: *Salmonella* spp.

Note: For safety and biohazard guidelines, refer to the “Safety” appendix in the PrepSEQ™ Rapid Spin Sample Preparation Kit *Salmonella* spp. (PN 4412848).

Product overview

The PrepSEQ™ Rapid Spin Sample Preparation Kit provides a simple way to prepare DNA from *broth* cultures.

Kit contents

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

Table 1

Item	Quantity or volume
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

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

Storage

- Lysis Buffer: 4 °C
- Spin columns and tubes: 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.

Overview

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

- *Salmonella* spp. enrichment
- Sample preparation

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

For some foods with a high lipid content, such as baby formula, soft cheese, whole milk, smoked salmon (lox), and chicken wing samples, the PrepSEQ™ Rapid Spin Extra Clean protocol is recommended.

For information on preparing:

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

Before you begin

Before starting your sample extraction:

- Set the block heater temperature to 95 °C.
- Label 1.5-mL microcentrifuge tubes.

Kit workflow

The figure below shows a sample processing workflow.

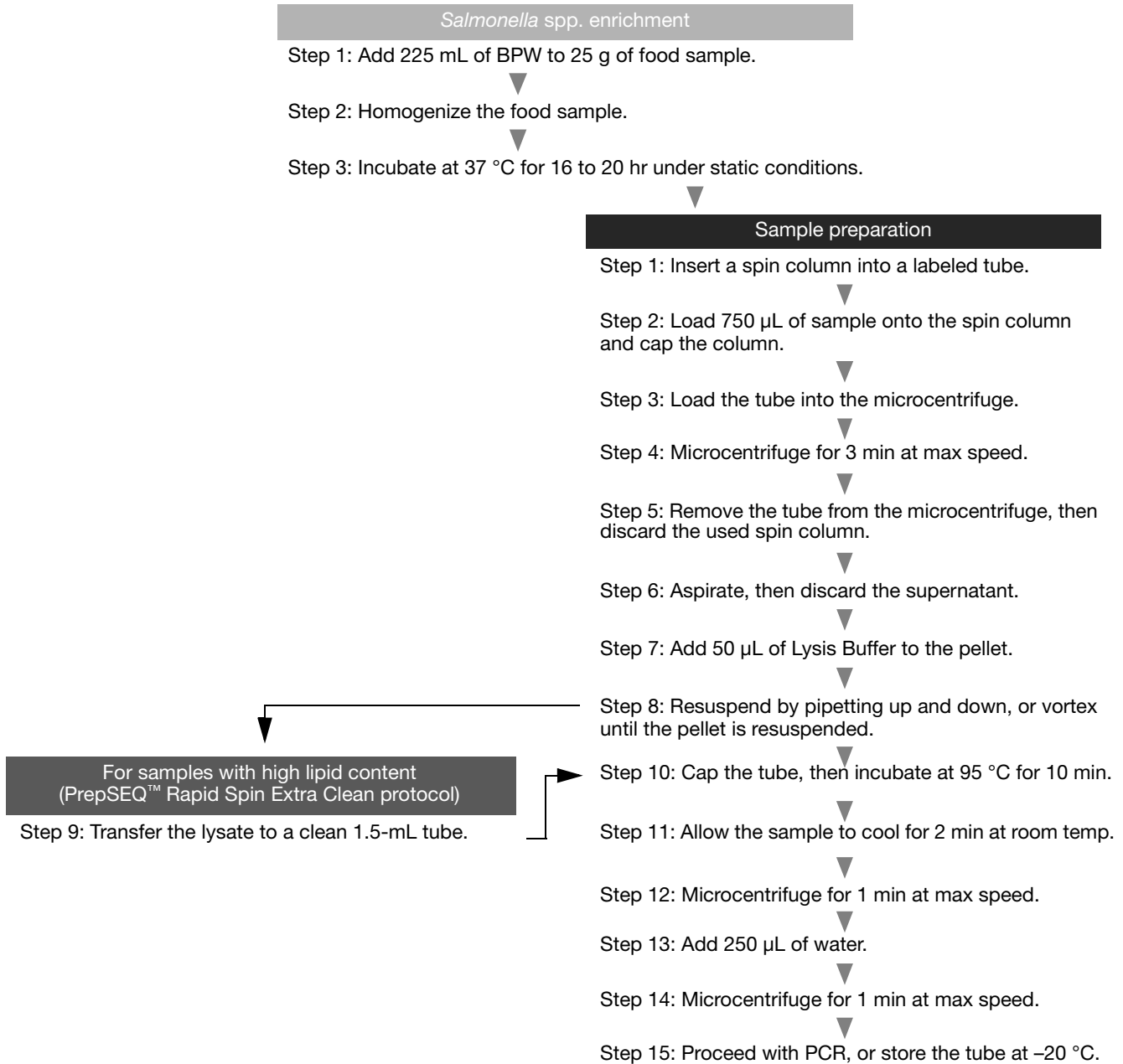


Figure 1 Food sample preparation using the PrepSEQ™ Rapid Spin Sample Preparation Kit

Notes

© Copyright 2010, Life Technologies Corporation. All rights reserved. **For Research Use Only. Not for use in diagnostic procedures.**

NOTICE TO PURCHASER: PLEASE REFER TO THE PREPSEQ NUCLEIC ACID EXTRACTION KIT PRODUCT INSERT AND PROTOCOL FOR LIMITED LABEL LICENSE OR DISCLAIMER INFORMATION.

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

05/2010

Part Number 4412849 Rev. B 05/2010



Headquarters

5791 Van Allen Way | Carlsbad, CA 92008 USA
Phone 760.603.7200
www.lifetechnologies.com

Technical Resources and Support

For the latest technical resources and support information for all locations, please refer to our Web site at www.appliedbiosystems.com/support