Ion Plus Fragment Library Kit

Pub. Part no. 4471992A Rev. Date 14 September 2011

Description

The Ion Plus Fragment Library Kit contains one box:

• Ion Plus Fragment Library Kit Part no. 4471252

The Ion Plus Fragment Library Kit contains sample preparation reagents to prepare up to 10 DNA libraries for semiconductor-based sequencing. Store the kit components as directed in the table. Please refer to the *Ion Xpress*TM *Plus gDNA and Amplicon Library Preparation User Guide* (Part no. 4471989) for more information.

IMPORTANT! Do not substitute components in the Ion Plus Fragment Library Kit with components in the Ion Fragment Library Kit (Part no. 4466464).

Precautions

Handle biological specimens and all other materials according to local laboratory guidelines and regulations. Avoid crosscontamination of samples and reagents to ensure quality results. If equipment or reagents are used in a manner not specified by the manufacturer, the protection provided by the equipment may be impaired.

For safety and biohazard guidelines, refer to the "Safety" section in the *Ion Xpress*[™] *Plus gDNA and Amplicon Library Preparation User Guide*. For every chemical, read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Note: Refer to the expiration date on the package label. Do not use kits after the expiration date.

Contact information

Phone: In North America: 1-87-SEQUENCE (1-877-378-3623)

Outside of North America: +1-203-458-8552

Address:

246 Goose Lane	7000 Shoreline Court	
Suite 100	Suite 201	
Guilford, CT 06437	South San Francisco, CA 94080	
USA	USA	

Web site: www.iontorrent.com

Ion community: ioncommunity.iontorrent.com Service email: ionsupport@lifetech.com

Kit components

Ion Plus Fragment Library Kit (Part no. 4471252)				
Components	Cap color	Quantity	Volume	Storage
5× End Repair Buffer	Red	1 tube	400 µL	
End Repair Enzyme	Orange	1 tube	20 µL	
10× Ligase Buffer	Yellow	1 tube	200 µL	
DNA Ligase	Blue	1 tube	40 µL	
Adapters	Green	1 tube	100 µL	
Platinum® PCR SuperMix High Fidelity	Black	2 tubes	1 mL	-30°C to -10°C
Library Amplification Primer Mix	White	1 tube	100 µL	
Nick Repair Polymerase	Clear	1 tube	160 µL	
dNTP Mix	Violet	1 tube	40 µL	
Low TE	Clear	2 tubes	1.5 mL	15 to 30°C or -30°C to -10°C



For Research Use Only. Not intended for any animal or human therapeutic or diagnostic use. COPYRIGHT

© 2011, Life Technologies Corporation. All rights reserved. No part of this publication may be reproduced, transmitted, transcribed, stored in retrieval systems, or translated into any language or computer language, in any form or by any means: electronic, mechanical, magnetic, optical, chemical, manual, or otherwise, without prior written permission from Life Technologies Corporation (hereinafter, "Life Technologies"). The information in this guide is subject to change without notice. Life Technologies and/or its affiliates reserve the right to change products and services at any time to incorporate the latest technological developments. Although this guide has been prepared with every precaution to ensure accuracy, Life Technologies and/or its affiliates assume no liability for any errors or omissions, nor for any damages resulting from the application or use of this information. Life Technologies welcomes customer input on corrections and suggestions for improvement.

TRADEMARKS

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

LIMITED LICENSE

All rights in the Personal Genome MachineTM (PGMTM) and in the information provided herein are owned by Life Technologies and/or its affiliates. Subject to the terms and conditions that govern your use of such products and information, Life Technologies and/or its affiliates grant you a nonexclusive, nontransferable, non-sublicensable license to use such products and information only in accordance with the manuals and written instructions provided by Life Technologies and/or its affiliates. You understand and agree that except as expressly set forth in the terms and conditions governing your use of such products, that no right or license to any patent or other intellectual property owned or licensable by Life Technologies and/or its affiliates is conveyed or implied by providing these products. In particular, no right or license is conveyed or implied to use these products in combination with any product not provided or licensed to you by Life Technologies and/or its affiliates for such use.

Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

