

Ion Shear™ Plus Reagents Kit

Pub. Part no. 100013885

MAN0005201

Rev. Date 22 November 2011

Kit components

The Ion Shear™ Plus Reagents Kit (Part no. 4471248) contains one box:

Ion Shear™ Plus Reagents Kit (Part no. 4471248)				
Components	Cap color	Quantity	Volume	Storage
Ion Shear™ Plus Enzyme Mix II	Clear	2 tubes	2 × 100 µL	-30°C to -10°C
Ion Shear™ Plus 10X Reaction Buffer	Clear	2 tubes	2 × 50 µL	
Ion Shear™ Plus Stop Buffer	Clear	2 tubes	2 × 50 µL	

Description

The Ion Shear™ Plus Reagents Kit is designed for the enzymatic shearing of DNA for use in Ion library preparation workflows. This kit is a component of the Ion Xpress™ Plus Fragment Library Kit (Part no. 4471269). Refer to the *Ion Xpress™ Plus gDNA and Amplicon Library Preparation User Guide* (Part no. 4471989) for a detailed user protocol.

IMPORTANT! Do not substitute components in the Ion Shear™ Plus Reagents Kit with components in the Ion Shear™ Reagents Kit (Part no. 4468655).

Precautions

Handle biological specimens and all other materials according to local laboratory guidelines and regulations. Avoid cross-contamination 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
Suite 100
Guilford, CT 06437
USA

7000 Shoreline Court
Suite 201
South San Francisco, CA 94080
USA

Web site: www.iontorrent.com

Ion community: ioncommunity.iontorrent.com

Support email: ionsupport@lifetech.com

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 Machine™ (PGM™) 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, non-transferable, 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.

**For support visit www.iontorrent.com/support
or email ionsupport@lifetech.com**

www.lifetechnologies.com

