


Vortex Adapter for Vortex Genie® 2 Mixer

Catalog Number AM10024

Pub. No. 4386453 Rev. B

| Contents | Quantity |
|--|----------|
| Vortex Adapter for Vortex Genie® 2 Mixer | 1 |

 **WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from www.lifetechnologies.com/support.

Product description

The Vortex Adapter attaches to Vortex Genie® 2 (Scientific Industries) vortex mixers and holds up to 12 microcentrifuge tubes for hands-free agitation.

Using the Vortex Adapter

To install on vortex mixers, align the flat surface inside the black adapter on the bottom with the flat surface on the vortex mixer head, and press the Vortex Adapter down until it snaps on. To remove the Vortex Adapter, place thumb on the label, with fingers underneath platform, then pull up and away.

Applications: The Vortex Adapter is required for use with the RiboPure™-Bacteria Kit (Cat. no. AM1925). During the RiboPure™-Bacteria procedure, position the sample tubes horizontally on the vortex adapter by sliding the tubes with the tube caps towards the center until they click; always oppose for balance. Vortex at maximum speed for 10 min.

For additional information on the procedure and protocol, please refer to the RiboPure™-Bacteria Kit Instruction Manual, found on the web at www.lifetechnologies.com.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

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

The information in this guide is subject to change without notice.

DISCLAIMER: LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) 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 USE LABEL LICENSE: Research Use Only : Notice to Purchaser: The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of 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.

© 2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners. Vortex Genie is a registered trademark of Scientific Industries, Inc.

For support visit lifetechnologies.com/support or email techsupport@lifetech.com

lifetechnologies.com