

## Preparing Small RNA Libraries for PGM™ Sequencing

Publication No. MAN0006951

Rev. Date: 28 June 2012

### RNA Sequencing

The Total Exosome RNA and Protein Isolation Kit (Cat. no. 4478545) is recommended for purification of exosomal RNA. RNA content of exosomes derived from at least 5 mL of cell media using the Total Exosome Isolation (from cell culture media) reagent or at least 200 µL of serum with the Total Exosome Isolation (from serum) reagent can be sequenced using the techniques described below.

For exosomes isolated using ultracentrifugation protocols (with or without sucrose gradient), we recommend increasing the input volume by 10-fold, because of very low recovery with this method.

For the complete protocol, refer to the manual for the Total Exosome RNA and Protein Isolation Kit at [www.lifetechnologies.com](http://www.lifetechnologies.com).

Small RNA libraries can be conveniently prepared with the Ion Total RNA-Seq Kit v2 (Cat. no. 4475936) and sequenced on the Ion Torrent™ PGM™.

The following version of the Total RNA-Seq Kit protocol has been modified to accommodate these specific characteristics of exosome samples:

- Relatively low amount of RNA
- Majority of the RNA cargo is short <200 nt

**Note:** This protocol is just a guide, and further modifications may be required, depending on the objectives of your study, the origin of the exosomes, the concentration and purity of RNA, the availability of certain kits/reagents, and other factors. For additional details on small RNA library construction, refer to the Ion Total RNA-Seq Kit v2 User Guide at [www.lifetechnologies.com](http://www.lifetechnologies.com).

### Experimental Overview

#### (Optional) RNase III fragmentation

RNase III fragmentation is recommended when performing analysis of long RNA. However, it is not necessary if the primary goal is analysis of short RNA (<200 nt). For additional details on fragmentation clean up, refer to the manual for the Purelink® RNA Micro Kit at [www.lifetechnologies.com](http://www.lifetechnologies.com).

1. Incubate the reaction in a thermal cycler at 37°C for 30 seconds.
2. Clean up 100 µL of the fragmented RNA with the Purelink® RNA Micro Kit (Cat no. 12183-016).
3. Add 100 µL of Lysis Buffer and 400 µL of 100% ethanol, and then mix well.
4. Bind, wash, and elute the RNA (in 12 µL) from the column.

### Small Library Construction

#### Hybridize and ligate the small RNA

1. Hybridize at 65°C for 10 minutes and 16°C for 5 minutes.
2. Perform ligation by incubating at 16°C for 16 hours (overnight).

#### Perform reverse transcription

1. Incubate the reverse transcription (RT) mix with the ligated RNA sample at 70°C for 10 minutes.
2. Perform RT reaction at 42°C for 30 minutes.

#### Purify cDNA using MagMAX™ Beads

1. Add 5 µL beads to one well of a 96-well plate for each sample.
2. Add 250 µL Binding Solution Concentrate to each well containing beads, and pipet up and down 10 times.
3. Add 60 µL of nuclease-free water to each of the 40 µL RT reactions and transfer to one of the wells of the 96-well plate.
4. Add 275 µL 100% ethanol to each well and mix by pipetting up and down.
5. Wash and elute cDNA (in 12 µL) from MagMAX™ beads.

#### Amplify the cDNA

1. Perform 18 cycles of PCR (total) with Platinum® PCR SuperMix High Fidelity.

#### Purify the amplified DNA

1. Add 5 µL beads to one well on a 96-well plate for each sample.
2. Add 280 µL Binding Solution Concentrate to each well containing beads, and pipet up and down 10 times.
3. Add 27 µL of nuclease-free water to each of the 53 µL PCR reactions.
4. Add 230 µL 100% ethanol to each well and mix by pipetting up and down.
5. Wash and elute cDNA (in 10 µL) in from MagMAX™ beads.

#### Analyze the cDNA

1. Run 1 µL of the sample on Agilent DNA High Sensitivity chip to assess the yield and size distribution.
2. Determine the molar concentration of the library with the Agilent® 2100 Bioanalyzer® Instrument Expert software.

#### Perform sequencing on the Ion PGM™ System

1. Enter settings and run sample. Small RNA libraries should run at 160 flows (40 cycles).
2. Enter the remaining information as needed, then follow the remaining prompts to start the run.

## Limited Use Label License: Research Use Only

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](mailto:outlicensing@lifetech.com) or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

## 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](http://www.lifetechnologies.com/termsandconditions). If you have any questions, please contact Life Technologies at [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support).

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.

---

©2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

For support visit [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support) or email [techsupport@lifetech.com](mailto:techsupport@lifetech.com)



[www.lifetechnologies.com](http://www.lifetechnologies.com)