

Preparation and Sequencing of RNA Libraries with the Ion Personal Genome Machine® (PGM™) System

Publication Part Number 4478119 (Rev. C) Revision Date 17 May 2012

■ Quantitate the library by Bioanalyzer™ Instrument analysis	1
Required materials	1
Procedure	2
■ Determine the Template Dilution Factor of the RNA library	4
Determine the Template Dilution Factor to prepare template-positive ISPs on the Ion OneTouch™ System for whole transcriptome and small RNA libraries	4
Determine the Template Dilution Factor to prepare template-positive ISPs with the IKA® Ultra-Turrax® Tube Drive	4
■ Set up the Ion PGM™ System for RNA libraries	5

Quantitate the library by Bioanalyzer™ Instrument analysis

Quantitate the library to determine its molar concentration. After determining the molar concentration of the library, calculate the Template Dilution Factor to dilute the library for preparation of template-positive Ion Sphere™ Particles (see [“Determine the Template Dilution Factor of the RNA library” on page 4](#)).

Note: For complete procedures including safety on RNA library preparation, refer to the *Ion Total RNA-Seq Kit v2 User Guide* (Pub. no. 4476286).

Required materials **Note:** For all required materials, consumables, and equipment, refer to the appropriate user guide.

Required kits

Item†	Source/Cat. no.
Agilent® DNA 1000 Kit	Agilent 5067-1504
Agilent® High Sensitivity DNA Kit	Agilent 5067-4626

† For the Safety Data Sheet (SDS) of any chemical not distributed by Life Technologies, contact the chemical manufacturer. Before handling any chemicals, refer to the SDS provided by the manufacturer, and observe all relevant precautions.

Optional kits

Item†	Source/Cat. no.
Qubit® dsDNA HS Assay Kit	Life Technologies Q32851

† For the Safety Data Sheet (SDS) of any chemical not distributed by Life Technologies, contact the chemical manufacturer. Before handling any chemicals, refer to the SDS provided by the manufacturer, and observe all relevant precautions.

Required equipment

Item	Source/Cat. no.
Agilent® 2100 Bioanalyzer™ instrument	Agilent G2938A
NanoDrop® Spectrophotometer	Thermo Scientific

Optional equipment

Item	Source/Cat. no.
Qubit® 2.0 Fluorometer	Life Technologies Q32866

Required consumables

Item	Source
(Optional) Low TE Buffer	Major Laboratory Supplier (MLS)

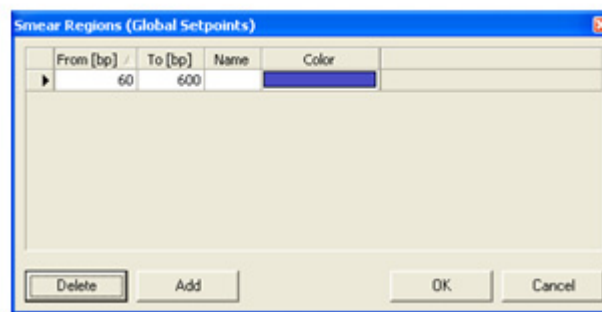
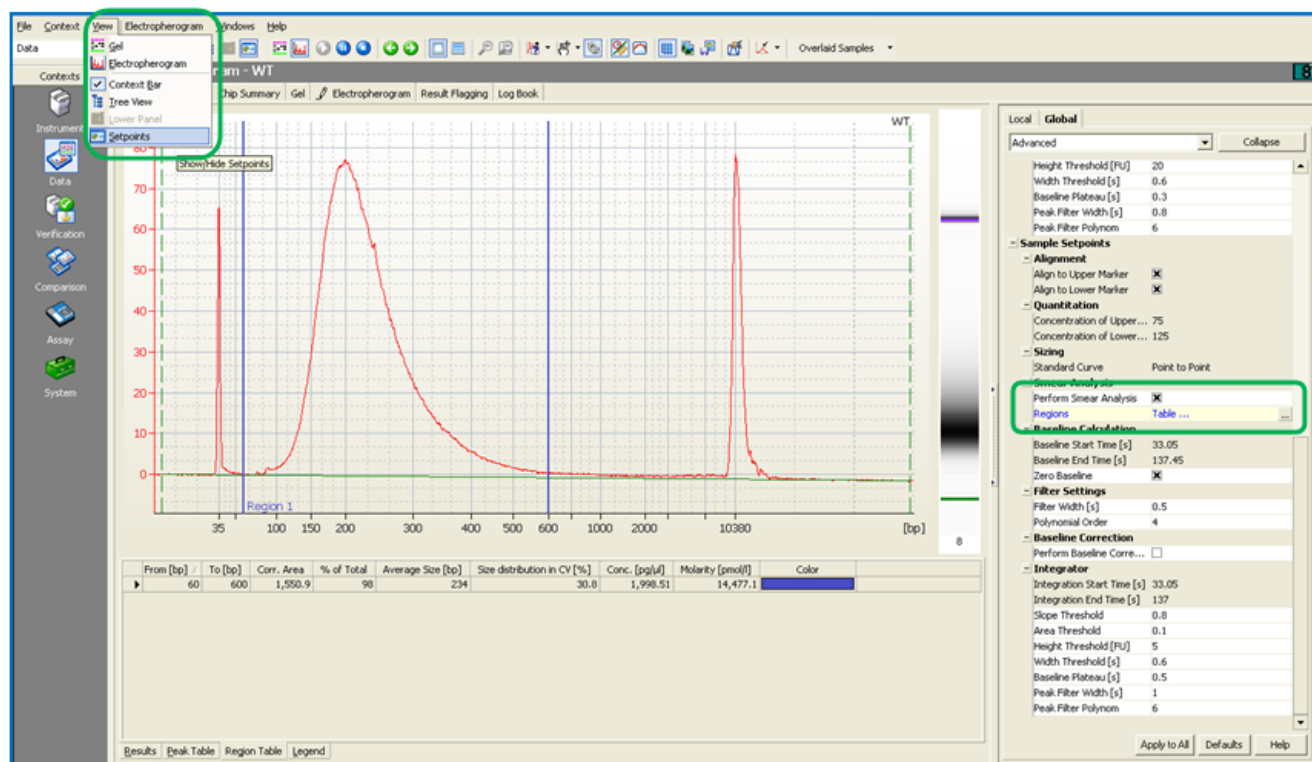
Procedure

1. Quantitate the purified library using 1 µL of sample with a Qubit® 2.0 Fluorometer, NanoDrop® Spectrophotometer, or other UV spectrophotometer (OD 260 nm).
2. Analyze 1 µL of the library using the appropriate chip on the Agilent® 2100 Bioanalyzer™ Instrument. If the library concentration is:
 - 1–50 ng/µL: Use the Agilent® DNA 1000 Kit.
 - 5–1000 pg/µL: Use the Agilent® High Sensitivity DNA Kit.
3. Determine the molar concentration of the library with the Agilent® 2100 Bioanalyzer™ Instrument Expert software: Perform a smear analysis by following the manufacturer's instructions. In the Bioanalyzer™ Instrument Expert software, select: **View ▶ Setpoints ▶ Smear Analysis**.

Example

The following figure shows a 60–600 bp region set for smear analysis. The software automatically calculates the mass concentration, either in:

- ng/ μ L and displays nMol/L with the Agilent® DNA 1000 Kit
or
- pg/ μ L and displays pmol/L with the Agilent® High Sensitivity DNA Kit:



Determine the Template Dilution Factor of the RNA library

The Template Dilution Factor is calculated from the molar concentration from Bioanalyzer™ Instrument analysis (see “[Quantitate the library by Bioanalyzer™ Instrument analysis](#)” on page 1). Calculate the Template Dilution Factor according to method. Next, use the Ion OneTouch™ 200 Template Kit v2 or the Ion PGM™ 200 Xpress™ Template Kit to prepare template-positive Ion Sphere™ Particles.

Note: For complete procedures on preparation of template-positive Ion Sphere™ Particles (ISPs) by the Ion OneTouch™ System, refer to the *Ion OneTouch™ 200 Template Kit v2 User Guide for use with the Ion OneTouch™ System* (Pub. no. 4478372).

Determine the Template Dilution Factor to prepare template-positive ISPs on the Ion OneTouch™ System for whole transcriptome and small RNA libraries

For the Template Preparation procedure using the Ion OneTouch™ 200 Template Kit v2, determine the Template Dilution Factor that gives 210×10^6 molecules of template per 20 μL for both whole transcriptome and small RNA libraries. Use a conversion factor of $8.3 \text{ nM} = 5 \times 10^9 \text{ molecules}/\mu\text{L}$, and the appropriate formula:

Template Dilution Factor =

$$(\text{Library concentration in nM}) \times [(5 \times 10^9 \text{ molecules}/\mu\text{L}) / (8.3 \text{ nM})] \times [20 \mu\text{L} / (210 \times 10^6 \text{ molecules})]$$

If your library concentration is $>20 \text{ nM}$, serially dilute 1:10 or 1:100 until your library concentration is $\leq 20 \text{ nM}$. (This step helps prevent a Template Dilution Factor of several thousands).

Example: If the library concentration is 10 nM :

Template Dilution Factor =

$$(10 \text{ nM}) \times [(5 \times 10^9 \text{ molecules}/\mu\text{L}) / (8.3 \text{ nM})] \times [20 \mu\text{L} / (210 \times 10^6 \text{ molecules})] = 574$$

Thus, 1 μL of library mixed with 573 μL of Nuclease-Free Water (1:574 dilution) yields approximately 210×10^6 molecules per 20 μL .

Determine the Template Dilution Factor to prepare template-positive ISPs with the IKA® Ultra-Turrax® Tube Drive

Note: For complete procedures on preparation of template-positive Ion Sphere™ Particles (ISPs) with the IKA® Ultra-Turrax® Tube Drive, refer to the *Ion PGM™ 200 Xpress™ Template Kit User Guide* (Pub. no. 4474291).

For the Template Preparation procedure using the Ion PGM™ 200 Xpress™ Kit, determine the Template Dilution Factor that gives 315×10^6 molecules of template per 18 μL for both whole transcriptome and small RNA libraries. Use a conversion factor of $8.3 \text{ nM} = 5 \times 10^9 \text{ molecules}/\mu\text{L}$, and the appropriate formula:

Template Dilution Factor =

$$(\text{Library concentration in nM}) \times [(5 \times 10^9 \text{ molecules}/\mu\text{L}) / (8.3 \text{ nM})] \times [18 \mu\text{L} / (315 \times 10^6 \text{ molecules})]$$

If your library concentration is $>20 \text{ nM}$, serially dilute 1:10 or 1:100 until your library concentration is less than or equal to 20 nM . (This step helps prevent a Template Dilution Factor of several thousands).

Example: The RNA library concentration is 400 nM

Dilute 1 μL of library in 99 μL of Nuclease-Free Water (1:100 dilution) to yield a final library concentration of 4 nM . Use this final library concentration to calculate the Template Dilution Factor:

Template Dilution Factor =
 $(4 \text{ nM}) \times [(5 \times 10^9 \text{ molecules}/\mu\text{L}) / (8.3 \text{ nM})] \times [18 \mu\text{L} / (315 \times 10^6 \text{ molecules})] = 138$

Thus, 1 μL of the 4-nM library dilution mixed with 137 μL of Nuclease-Free Water (1:138 dilution) yields approximately 315×10^6 molecules per 18 μL . Use 18 μL of the final library dilution to prepare the aqueous master mix.

Set up the Ion PGM™ System for RNA libraries

1. Enter these settings on the Run Info screen of the Ion PGM™ Sequencer:

If the touchscreen prompt says...	Select...
Runtype	GENS
Library	Appropriate library reference
Barcode	• <i>Non-barcoded</i> libraries: RNA_Barcode_None • <i>Barcoded</i> libraries prepared with the Ion Xpress™ RNA-Seq Barcode 01–16 Kit: IonXpressRNA
Number of flows	• <i>Whole transcriptome</i> libraries: 520 flows (130 cycles) • <i>Small RNA</i> libraries: 160 flows (40 cycles) Note: To change cycling conditions for different read lengths, enter a different number of flows.

IMPORTANT! To insure proper trimming of the sequences prior to alignment with a reference, for all libraries prepared with the Ion Total RNA-seq Kit v2, you must select an option from the barcode pull-down menu. If the RNA-seq library does *not* have barcodes, select **RNA_Barcode_None** in the barcode dialog box. If the library was prepared with the Ion Xpress™ RNA-Seq Barcode 01–16 Kit, select **IonXpressRNA**.

2. Enter the remaining information as needed.
3. Click **Next**, then follow the remaining prompts to start the run.

For research use only. Not intended for diagnostic procedures.

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

DISCLAIMERS

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 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, 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 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 or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

TRADEMARKS

The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliates or their respective owners. Bioanalyzer is a trademark of Agilent Technologies, Inc. Agilent is a registered trademark of Agilent Technologies, Inc. IKA and Ultra-Turrax are registered trademarks of IKA Werke GmbH & Co. NanoDrop is a registered trademark of NanoDrop Technologies LLC. Microsoft and Excel are registered trademarks of Microsoft Corporation.

© 2012, Life Technologies Corporation. All rights reserved.

Headquarters

5791 Van Allen Way | Carlsbad, CA 92008 USA

Phone +1 760 603 7200 | Toll Free in USA 800 955 6288

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

www.lifetechnologies.com

