

SureSelect^{XT} Automated RNA Target Enrichment

**Automated Poly-A Selection,
Strand-Specific mRNA Library
Preparation and Target Enrichment for
the Illumina Platform**

Protocol

Version G0, June 2020

**SureSelect platform manufactured with Agilent
SurePrint Technology**

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



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2013, 2014, 2017, 2018, 2020

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

Manual Part Number

G9691-90020

Edition

Version G0, June 2020

Printed in USA

Agilent Technologies, Inc.
5301 Stevens Creek Blvd
Santa Clara, CA 95051 USA

Acknowledgment

Oligonucleotide sequences © 2006, 2008, and 2011 Illumina, Inc. All rights reserved. Only for use with the Illumina sequencer systems and associated assays.

Technical Support

For US and Canada

Call (800) 227-9770 (option 3,4,4)

Or send an e-mail to:

ngs.support@agilent.com

For all other regions

Agilent's world-wide Sales and Support Center contact details for your location can be obtained at

www.agilent.com/en/contact-us/page.

Warranty

The material contained in this document is provided "as is," and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

Restricted Rights Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

Notice to Purchaser: Limited License

Limited Use Label License: This product and its use are the subject of one or more issued and/or pending U.S. and foreign patent applications owned by Max Planck Gesellschaft exclusively licensed to New England Biolabs, Inc. and sublicensed to Agilent Technologies. The purchase of this product from Agilent Technologies, Inc., its affiliates, or its authorized resellers and distributors conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for profit entity). The purchase of this product does not convey a license under any claims in the foregoing patents or patent applications directed to producing the product. The buyer cannot sell or otherwise transfer this product or its components to a third party or otherwise use this product for the following COMMERCIAL PURPOSES: (1) use of the product or its components in manufacturing; or (2) use of the product or its components for therapeutic or prophylactic purposes in humans or animals.

Safety Notices

CAUTION

A **CAUTION** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a **CAUTION** notice until the indicated conditions are fully understood and met.

WARNING

A **WARNING** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

In this Guide...

This guide describes an optimized protocol to prepare target-enriched mRNA sequencing libraries from total RNA samples using the SureSelect Automated Strand-Specific RNA Library Prep and Target Enrichment system.

This protocol is specifically developed for RNA library preparation for Illumina paired-end multiplexed sequencing. Sample processing steps are automated using the NGS Workstation.

1 Before You Begin

This chapter contains information (such as procedural notes, safety information, required reagents and equipment) that you should read and understand before you start an experiment.

2 Using the Agilent NGS Workstation for SureSelect RNA Library Preparation

This chapter contains an orientation to the Agilent NGS Workstation, an overview of the SureSelect RNA protocol, and considerations for designing SureSelect experiments for automated processing using the Agilent NGS Workstation.

3 Sample Preparation

This chapter describes the steps to prepare strand-specific mRNA libraries from total RNA samples for sequencing on the Illumina platform.

4 Hybridization

This chapter describes the steps to hybridize the prepped library with the SureSelect or ClearSeq RNA Capture Library and to capture the targeted sequences.

5 Indexing and Sample Prep for Multiplexed Sequencing

This chapter describes the steps to index by amplification, purify, and assess the quality and quantity of the target-enriched samples. Indexed samples are pooled by mass prior to sequencing.

6 Reference

This chapter contains reference information.

What's New in Version G0

- Support for renamed Library Preparation Kit components. See [Table 75](#) on page 134 and [Table 76](#) on page 135 for a summary of component name changes. The new component names are utilized throughout the protocols in this document. No changes were made to the reagent formulations or methods of use.
- Support for revised Poly-A Selection Module subkit. See [Table 75](#) on page 134 and [Table 77](#) on page 136 for a summary of component name and configuration changes. The new component names are utilized throughout the protocols in this document. No changes were made to the reagent formulations or methods of use. Water is no longer provided in this module (see [Table 77](#) on page 136) and has been added to the list of required reagents in [Table 1](#) on page 12.
- Updates to ordering information for custom RNA target enrichment probes (see [Table 2](#) on page 13) and updates to probe nomenclature in the Hybridization protocol steps (see [page 87](#) to [page 90](#)).
- Updates to sequencing support guidelines for RNA strandedness (see [page 131](#)).
- Updated recommendation for control reference RNA (see [Table 4](#) on page 16).
- Updates to ordering information for 1X TE Buffer and AMPure XP Kits (see [Table 1](#) on page 12).
- Updated Technical Support contact information (see [page 2](#)).
- Updated *Notice to Purchaser* (see [page 2](#)).
- End of support for component kit 5190-4394. This kit was replaced with component kit 5190-8646 in May 2018 for use in the current on-bead post-capture PCR method. See [Table 75](#) on page 134 and [Table 78](#) on page 136 for current kit contents.

Content

1	Before You Begin	9
	Procedural Notes	10
	Safety Notes	11
	Required Reagents	12
	Required Equipment	14
	Optional Reagents and Equipment	16
2	Using the Agilent NGS Workstation for SureSelect RNA Library Preparation	17
	About the Agilent NGS Workstation	18
	About the Bravo Platform	18
	VWorks Automation Control Software	22
	Overview of the SureSelect RNA Library Prep Procedure	28
	Experimental Setup Considerations for Automated Runs	30
	Considerations for Placement of RNA Samples in 96-well Plates for Automated Processing	31
	Considerations for Equipment Setup	31
	PCR Plate Type Considerations	32
3	Sample Preparation	33
	Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA	34
	Step 2. Purify first-strand cDNA using AMPure XP beads	55
	Step 3. Prepare cDNA libraries for Hybridization	59
	Step 4. Amplify cDNA libraries by PCR	67
	Step 5. Purify amplified DNA using AMPure XP beads	75
	Step 6. Assess library DNA quantity and quality	79

Contents

4	Hybridization	83
	Step 1. Aliquot prepped DNA libraries for hybridization	84
	Step 2. Hybridize the DNA library to the Probe	87
	Step 3. Capture the hybridized DNA	100
5	Indexing and Sample Prep for Multiplexed Sequencing	109
	Step 1. Amplify the captured libraries to add index tags	110
	Step 2. Purify the amplified indexed libraries using AMPure XP beads	119
	Step 3. Assess DNA quality and quantity	123
	Step 4. Pool samples for multiplexed sequencing	127
	Step 5. Prepare and analyze sequencing samples	129
6	Reference	133
	Kit Contents	134
	Nucleotide Sequences of SureSelect ^{XT} Indexes A01 to H12	138
	Plasticware quantities for automation protocols	139



1 Before You Begin

Procedural Notes	10
Safety Notes	11
Required Reagents	12
Required Equipment	14
Optional Reagents and Equipment	16

Make sure you read and understand the information in this chapter and have the necessary equipment and reagents listed before you start an experiment.

NOTE

This protocol describes automated sample processing using the Agilent NGS Workstation. For non-automated sample processing procedures for Agilent's SureSelect Strand-Specific RNA Target Enrichment Kit for sequencing on the Illumina platform, see publication G9691-90000.



Procedural Notes

- Certain protocol steps require the rapid transfer of sample plates between the Bravo deck and a thermal cycler. Locate your thermal cycler in close proximity to the Agilent NGS Workstation to allow rapid and efficient plate transfer.
- Use of Agilent's SureCycler 8800 thermal cycler and associated plasticware is recommended for optimal performance. The workflow is compatible with additional thermal cyclers, but performance should be validated before running a large number of samples. See [page 32](#) for a list of supported PCR plate types and ensure that the thermal cycler to be used is compatible with one of the supported PCR plate types.
- Prepare and load the Agilent NGS Workstation as detailed in each of the protocol steps before initiating each automated protocol run. When loading plates in the workstation's Labware MiniHub, always place plates in the orientation shown in [Figure 4](#) on page 39.
- To prevent contamination of reagents by nucleases, always wear powder-free laboratory gloves and use dedicated solutions and pipettors with nuclease-free aerosol-resistant tips.
- Avoid repeated freeze-thaw cycles of stock and diluted RNA and cDNA solutions. Possible stopping points, where samples may be stored at -20°C , are marked in the protocol. Do not subject the samples to multiple freeze/thaw cycles.
- When preparing master mix reagent stock solutions for use:
 - 1 Thaw the reagent vial as rapidly as possible without heating above room temperature.
 - 2 Mix thoroughly on a vortex mixer at high speed for 5 seconds, then briefly spin in a centrifuge to drive the contents off of walls and lid.
 - 3 Store vials used during an experiment on ice or in a cold block.
 - 4 Library Preparation Master Mixes should not be frozen and thawed more than five times. If you plan to use the reagents in more than five experiments, aliquot to multiple vials to minimize freeze/thaw cycles for each vial.
- In general, follow Biosafety Level 1 (BL1) safety rules.

Safety Notes

CAUTION

- Wear appropriate personal protective equipment (PPE) when working in the laboratory.
-

Required Reagents

Table 1 Required Reagents for SureSelect^{XT} RNA Target Enrichment Automation

Description	Vendor and part number
SureSelect or ClearSeq RNA Target Enrichment Probe	Select one probe library from Table 2
SureSelect ^{XT} RNA Reagent Kit*, Illumina platforms (ILM), 96 Samples*	Agilent p/n G9692B
Actinomycin D [†]	Sigma p/n A1410
DMSO	Sigma p/n D8418
Dynabeads M-270 Streptavidin Beads	Thermo Fisher Scientific p/n 65306
AMPure XP Kit	Beckman Coulter Genomics
5 mL	p/n A63880
60 mL	p/n A63881
450 mL	p/n A63882
1X Low TE Buffer (10 mM Tris-HCl, pH 7.5-8.0, 0.1 mM EDTA)	Thermo Fisher Scientific p/n 12090-015, or equivalent
100% Ethanol, molecular biology grade	Sigma-Aldrich p/n E7023
Nuclease-free Water (not DEPC-treated)	Thermo Fisher Scientific p/n AM9930

* Each 96-reaction kit contains sufficient reagents for 96 reactions used in runs that include at least 3 columns of samples per run.

† Actinomycin D should be obtained as a solid and prepared at 4 µg/µl concentration in DMSO then stored in single-use aliquots at –20°C, protected from light. The aliquots may be stored for up to one year before use. See [page 36](#) for more information.

CAUTION

Use only the recommended Dynabeads M-270 Streptavidin Beads for this automated protocol. Use of other streptavidin bead preparations may adversely affect performance and is not supported by Agilent.

Table 2 SureSelect and ClearSeq RNA Target Enrichment Probes

Capture Library	96 Samples
Custom RNA Target Enrichment Probes 1–499 kb	Please contact Agilent’s SureSelect support team or your local representative for assistance with custom RNA probe design and ordering.
Custom RNA Target Enrichment Probes 0.5–2.9 Mb	
Custom RNA Target Enrichment Probes 3– 5.9 Mb	
ClearSeq RNA Kinome XT	5190-7287

Required Equipment

Table 3 Required Equipment

Description	Vendor and part number
Agilent NGS Workstation Option B Contact Agilent Automation Solutions for more information: Customerservice.automation@agilent.com	Agilent p/n G5522A (VWorks software version 13.1.0.1366, 13.0.0.1360, or 11.3.0.1195) OR Agilent p/n G5574AA (VWorks software version 13.1.0.1366)
Bravo 96-well PCR plate insert (red)	Agilent p/n G5498B#13
Robotic Pipetting Tips (Sterile, Filtered, 250 µL)	Agilent p/n 19477-022
Thermal cycler and accessories	SureCycler 8800 Thermal Cycler (Agilent p/n G8810A), 96 well plate module (Agilent p/n G8810A) and compression mats (Agilent p/n 410187) or equivalent
PCR plates compatible with selected thermal cycler, e.g. Agilent semi-skirted PCR plate for the SureCycler 8800 Thermal Cycler When selecting plates for another thermal cycler, see page 32 for the list of PCR plates supported in automation protocols	Agilent p/n 401334
Eppendorf twin.tec full-skirted 96-well PCR plates	Eppendorf p/n 951020401 or 951020619
Thermo Scientific Reservoirs	Thermo Fisher Scientific p/n 1064-15-6
Nunc DeepWell Plates, sterile, 1.3-mL well volume	Thermo Fisher Scientific p/n 260251
Axygen 96 Deep Well Plate, 2 mL, Square Well (waste reservoirs; working volume 2.2 mL)	Axygen p/n P-2ML-SQ-C E & K Scientific p/n EK-2440
Nucleic acid surface decontamination wipes	DNA Away Surface Decontaminant Wipes, Thermo Fisher Scientific p/n 7008, or equivalent
Vacuum concentrator	Savant SpeedVac, model DNA120, with 96-well plate rotor, model RD2MP, or equivalent

Table 3 Required Equipment (continued)

Description	Vendor and part number
Magnetic separator	DynaMag-50 magnet, Thermo Fisher Scientific p/n 123-02D or equivalent
DNA Analysis Platform and Consumables	
Agilent 2100 Bioanalyzer Instrument	Agilent p/n G2939BA
Agilent 2100 Expert SW Laptop Bundle (optional)	Agilent p/n G2953CA
DNA 1000 Kit	Agilent p/n 5067-1504
High Sensitivity DNA Kit	Agilent p/n 5067-4626
OR	
Agilent 4200 TapeStation*	Agilent p/n G2991AA
96-well sample plates	Agilent p/n 5042-8502
96-well plate foil seals	Agilent p/n 5067-5154
8-well tube strips	Agilent p/n 401428
8-well tube strip caps	Agilent p/n 401425
D1000 ScreenTape	Agilent p/n 5067-5582
D1000 Reagents	Agilent p/n 5067-5583
High Sensitivity D1000 ScreenTape	Agilent p/n 5067-5584
High Sensitivity D1000 Reagents	Agilent p/n 5067-5585
Low-Adhesion Tubes (RNase, DNase, and DNA-free)	USA Scientific
1.5 mL	p/n 1415-2600
0.5 mL	p/n 1405-2600
P10, P20, P200 and P1000 pipettes	Pipetman P10, P20, P200, P1000 or equivalent
Ice bucket	
Powder-free gloves	
Vortex mixer	
Timer	

* DNA samples may also be analyzed using the 4150 TapeStation, p/n G2992AA. ScreenTape devices and associated reagents listed in this table are compatible with both platforms.

Optional Reagents and Equipment

Table 4 Optional Reagents and Equipment

Description	Vendor and part number
Labnet MPS1000 Mini Plate Spinner	Labnet International p/n C1000
Agilent QPCR Human Reference Total RNA	Agilent p/n 750500



2

Using the Agilent NGS Workstation for SureSelect RNA Library Preparation

About the Agilent NGS Workstation	18
Overview of the SureSelect RNA Library Prep Procedure	28
Experimental Setup Considerations for Automated Runs	30

This chapter contains an orientation to the Agilent NGS Workstation, an overview of the SureSelect strand-specific RNA library preparation and target enrichment protocol, and considerations for designing SureSelect RNA experiments for automated processing using the Agilent NGS Workstation.



About the Agilent NGS Workstation

CAUTION

Before you begin, make sure that you have read and understand operating, maintenance and safety instructions for using the Bravo platform and additional devices included with the workstation. Refer to the user guides listed in [Table 5](#).

Review the user guides listed in [Table 5](#) (available at Agilent.com) to become familiar with the general features and operation of the Agilent NGS Workstation Option B components. Instructions for using the Bravo platform and other workstation components for the SureSelect^{XT} RNA Target Enrichment workflow are detailed in this user guide.

Table 5 Agilent NGS Workstation components User Guide reference information

Device	User Guide part number
Bravo Platform	G5562-90000
VWorks Software	G5415-90068 (VWorks versions 13.1.0.1366 and 13.0.0.1360), or G5415-90063 (VWorks version 11.3.0.1195)
BenchCel Microplate Handler	G5400-90004
Labware MiniHub	G5471-90002
PlateLoc Thermal Microplate Sealer	G5402-90001

About the Bravo Platform

The Bravo platform is a versatile liquid handler with a nine plate-location platform deck, suitable for handling 96-well, 384-well, and 1536-well plates. The Bravo platform is controlled by the VWorks Automation Control software. Fitted with a choice of seven interchangeable fixed-tip or disposable-tip pipette heads, it accurately dispenses fluids from 0.1 µL to 250 µL.

Bravo Platform Deck

The protocols in the following sections include instructions for placing plates and reagent reservoirs on specific Bravo deck locations. Use [Figure 1](#) to familiarize yourself with the location numbering convention on the Bravo platform deck.

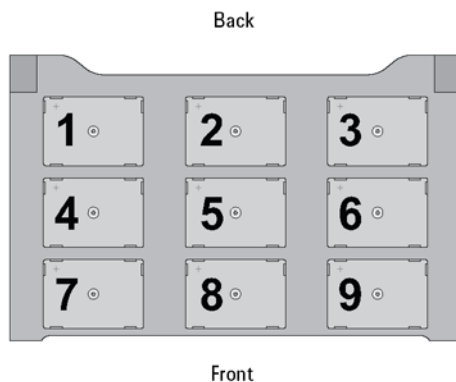


Figure 1 Bravo platform deck

Setting the Temperature of Bravo Deck Heat Blocks

Bravo deck positions 4 and 6 are equipped with Inheco heat blocks, used to incubate sample plates at defined temperatures during the run. Runs that include high- (85°C) or low- (4°C) temperature incubation steps may be expedited by pre-setting the temperature of the affected block before starting the run.

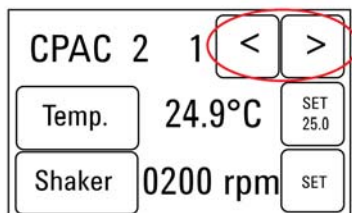
Bravo deck heat block temperatures may be changed using the Inheco Multi TEC Control device touchscreen as described in the steps below. See [Table 6](#) for designations of the heat block-containing Bravo deck positions on the Multi TEC control device.

Table 6 Inheco Multi TEC Control touchscreen designations

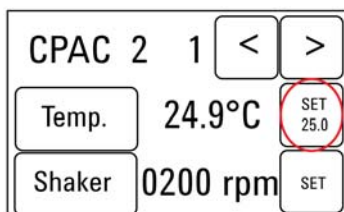
Bravo Deck Position	Designation on Inheco Multi TEC Control Screen
4	CPAC 2 1
6	CPAC 2 2

2 Using the Agilent NGS Workstation for SureSelect RNA Library Preparation About the Bravo Platform

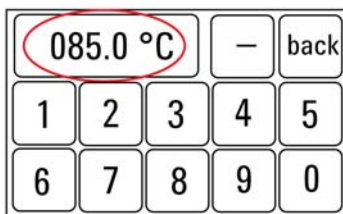
- 1 Using the arrow buttons, select the appropriate block (CPAC 2 block 1 or CPAC 2 block 2).



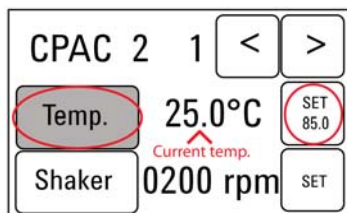
- 2 To set the temperature of the selected block, press the SET button.



- 3 Using the numeral pad, enter the desired temperature. The entered temperature appears in the top, left rectangle. Once the correct temperature is displayed, press the rectangle to enter the temperature.



- 4 Press the Temp button until the new temperature is displayed on the SET button and until the Temp button is darkened, indicating that the selected heat block is heating or cooling to the new temperature setting. The current temperature of the block is indicated in the center of the display.



Setting the Temperature of Bravo Deck Position 9 Using the ThermoCube Device

Bravo deck position 9 is equipped with a ThermoCube thermoelectric temperature control system, used to incubate components at a defined temperature during the run. During protocols that require temperature control at position 9, you will be instructed to start and set the temperature of the ThermoCube device before starting the run.

ThermoCube temperature settings are modified using the control panel (LCD display screen and four input buttons) on the front panel of the device using the following steps.

- 1 Turn on the ThermoCube and wait for the LCD screen to display **TEMP.**
- 2 Press the **UP** or **DOWN** button to change **SET TEMP 1** to the required set point.
- 3 Press the **START** button.

The ThermoCube will then initiate temperature control of Bravo deck position 9 at the displayed set point.

VWorks Automation Control Software

VWorks software, included with your Agilent NGS Workstation, allows you to control the robot and integrated devices using a PC. The Agilent NGS Workstation is preloaded with VWorks software containing all of the necessary SureSelect system liquid handling protocols. General instructions for starting up the VWorks software and the included protocols is provided below. Each time a specific VWorks protocol is used in the SureSelect procedure, any settings required for that protocol are included in the relevant section of this manual.

NOTE

The instructions in this manual are compatible with VWorks software version 13.1.0.1366, 13.0.0.1360 or 11.3.0.1195.

If you have questions about VWorks version compatibility, please contact service.automation@agilent.com.

Logging in to the VWorks software

- 1 Double-click the VWorks icon or the SureSelect_RNA_ILM.VWForm shortcut on the Windows desktop to start the VWorks software.
- 2 If User Authentication dialog is not visible, click **Log in** on the VWorks window toolbar.
- 3 In the User Authentication dialog, type your VWorks user name and password, and click **OK**. (If no user account is set up, contact the administrator.)

VWorks protocol and runset files

VWorks software uses two file types for automation runs, .pro (protocol) files and .rst (runset) files. Runset files are used for automated procedures in which the workstation uses more than one automation protocol during the run.

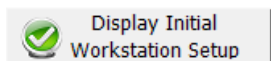
Using the SureSelect_RNA_ILM.VWForm to setup and start a run

Use the VWorks form SureSelect_RNA_ILM.VWForm, shown below, to set up and start each SureSelect automation protocol or runset.

The screenshot shows the 'SureSelect RNA for Illumina sequencers' VWorks form. It is divided into several sections:

- Parameters:**
 - 1) Select Protocol to Run: mRNA_Purification_v1.0.pro
 - AMPureXP Case: (empty)
 - 2) Select PCR Plate labware for Thermal Cycling: 96 ABI PCR half skirt in Red Alum Insert
 - 3) Select Number of Columns of Samples: 1
 - 4) Click button below to Display Initial Workstation Setup: Includes 'Display Initial Workstation Setup' and 'Clear Workstation Setup Display' buttons.
 - 5) Load labware according to Workstation Setup -->
- Controls:**
 - Once you have loaded labware according to Workstation Setup on right, click "Run Selected Protocol" to start run.
 - Buttons: Run Selected Protocol, Pause, Initialize all devices.
 - Full Screen, Gantt Chart, Elapsed Time: 00:00:00.
 - Reset All Form Selections to Defaults.
- Information:**
 - Currently Running Protocol: (empty)
- Workstation Setup:**
 - MiniHub:** A 5x4 grid table for MiniHub Cassette 1-4 and Shelf 1-5.
 - Bravo Deck:** A diagram of a Bravo deck with 9 positions: <Position 1>, <Position 2>, <Position 3>, <Pos 4: Peltier>, <Pos 5: Shaker>, <Pos 6: Peltier>, <Pos 7: Magnetic>, <Position 8>, <Pos 9: Chiller>.
 - BenchCel:** A 1x4 grid table for BenchCel Stacker 1-4.

- 1 Open the form using the SureSelect_RNA_ILM.VWForm shortcut on your desktop.
- 2 Use the drop-down menus on the form to select the appropriate SureSelect workflow step and number of columns of samples for the run.
- 3 Once all run parameters have been specified on the form, click **Display Initial Workstation Setup**.



2 Using the Agilent NGS Workstation for SureSelect RNA Library Preparation

VWorks Automation Control Software

- The Workstation Setup region of the form will then display the required placement of reaction components and labware in the NGS Workstation for the specified run parameters.

SureSelect RNA
for Illumina sequencers

Parameters

- Select Protocol to Run
mRNA_Purification_v1.0.pro
AMPureXP Case: Not Applicable
- Select PCR Plate labware for Thermal Cycling
96 ABI PCR half skirt in Red Alum Insert
- Select Number of Columns of Samples
1
- Click button below to Display Initial Workstation Setup
- Load labware according to Workstation Setup -->

Controls

Once you have loaded labware according to Workstation Setup on right, click "Run Selected Protocol" to start run.

Elapsed Time: 00:00:00

Information

Currently Running Protocol:

Workstation Setup

	MiniHub Cassette 1	MiniHub Cassette 2	MiniHub Cassette 3	MiniHub Cassette 4
Shelf 5	Empty Nunc DeepWell Plate	Empty Nunc DeepWell Plate	Empty Nunc DeepWell Plate	
Shelf 4		Bead Binding Buffer in twin.tec		
Shelf 3		Bead Elution Buffer in twin.tec		
Shelf 2	Empty Tip Box		Bead Wash Buffer in Nunc DeepWell	
Shelf 1	New Tip Box			Empty Tip Box

Bravo Deck

<Position 1> Waste Reservoir (Axygen 96DW)	<Position 2>	<Position 3>
<Pos 4: Peltier> _{RT} Oligo dT beads (Set labware in 2)	<Pos 5: Shaker>	<Pos 6: Peltier> _{RT} Empty Plate (Set labware in 2)
<Pos 7: Magnetic> Total RNA in twin.tec Plate	<Position 8>	<Pos 9: Chiller> _{0°C} Nunc Master Mix Plate (Col 1-2) on Silver Insert

BenchCel

BenchCel Stacker 1	BenchCel Stacker 2	BenchCel Stacker 3	BenchCel Stacker 4
1 Tip Box	Empty	Empty	Empty

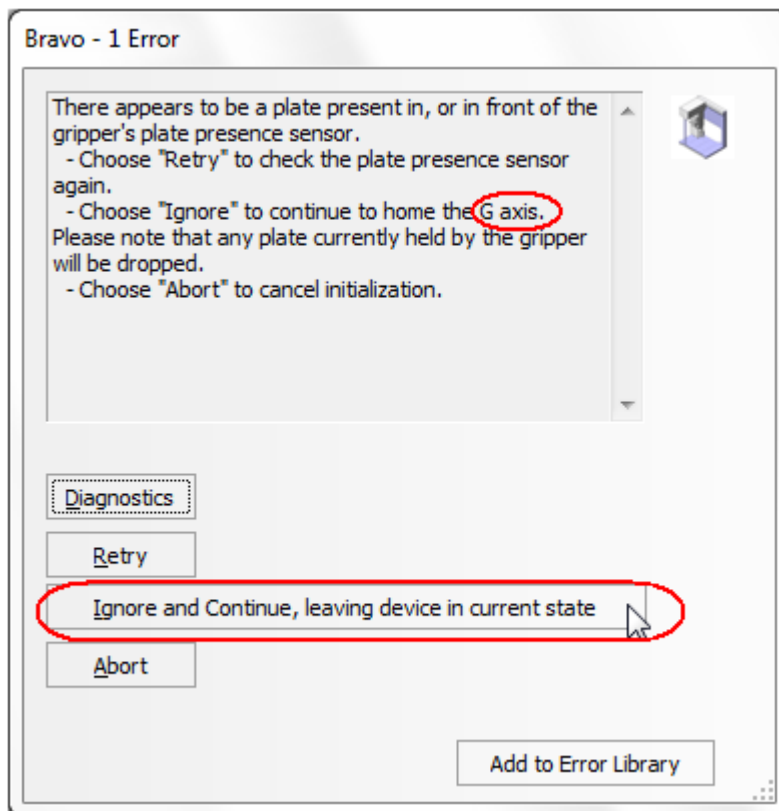
- After verifying that the NGS Workstation has been set up correctly, click **Run Selected Protocol**.



Error messages encountered at start of run

After starting the run, you may see the error messages displayed below. When encountered, make the indicated selections and proceed with the run. Encountering either or both of these error messages is not indicative of a problem with the NGS workstation or your run setup.

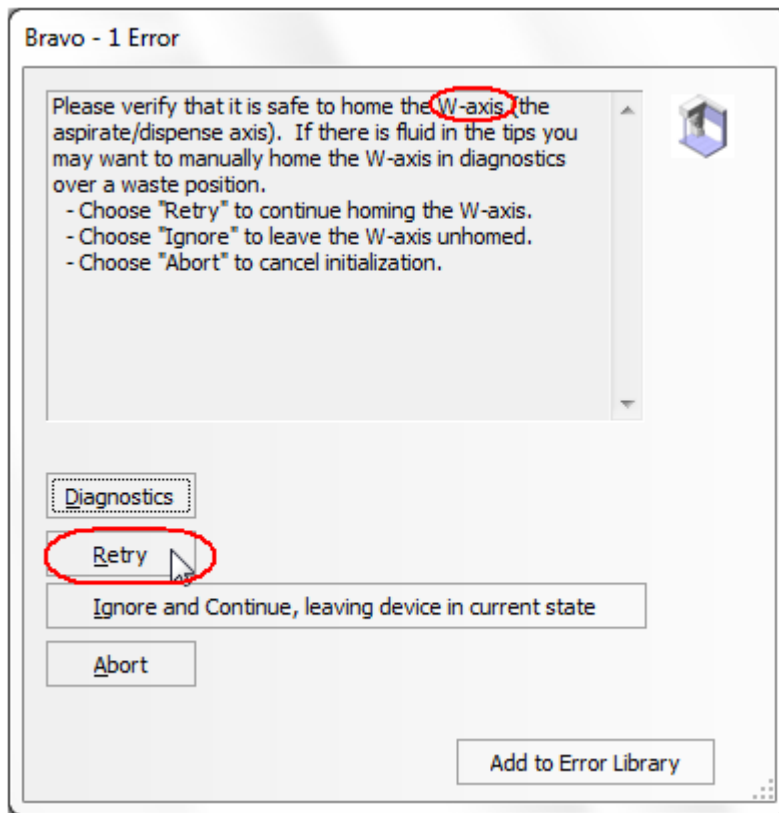
- 1 If you encounter the G-axis error message shown below, select **Ignore and Continue, leaving device in current state**.



2 Using the Agilent NGS Workstation for SureSelect RNA Library Preparation

VWorks Automation Control Software

- 2 If you encounter the W-axis error message shown below, select **Retry**.



Verifying the Simulation setting

VWorks software may be run in simulation mode, during which commands entered on screen are not completed by the NGS workstation. If workstation devices do not respond when you start a run, verify the simulation mode status in VWorks using the following steps.

- 1 Verify that **Simulation is off** is displayed on the status indicator (accessible by clicking **View > Control Toolbar**).



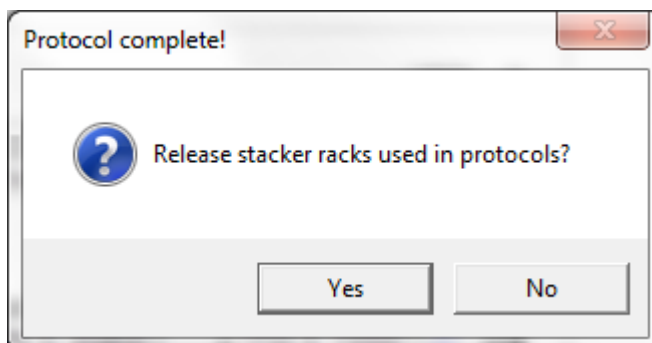
- 2 If the indicator displays **Simulation is on**, click the status indicator button to turn off the simulation mode.

NOTE

If you cannot see the toolbar above the SureSelect VWorks form, click the **Full Screen** button to exit full screen mode. If the toolbar is still not visible, right-click on the form and then select **Control Toolbar** from the menu.

Finishing a protocol or runset

The window below appears when each run is complete. Click **Yes** to release the BenchCel racks to allow removal of components used in the current run in preparation for the next .pro or .rst run.



Overview of the SureSelect RNA Library Prep Procedure

Figure 2 summarizes the SureSelect workflow for RNA samples to be sequenced using the Illumina sequencing platform. For each sample to be sequenced, an individual cDNA library is prepared. The libraries are then target enriched and tagged by PCR with an index sequence. Depending on the capacity of the sequencing platform, up to 96 samples can be pooled and sequenced in a single lane using the multiplex index tags that are provided with the SureSelect Strand-Specific RNA Library Prep kit.

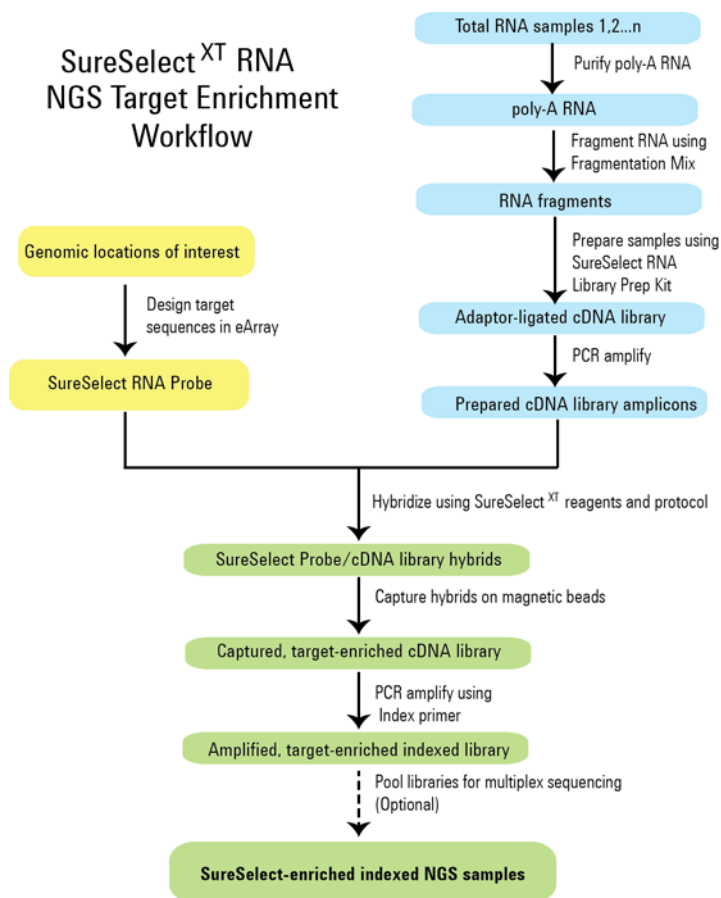


Figure 2 Overall sequencing sample preparation workflow

Table 7 summarizes how the VWorks protocols are integrated into the Strand-Specific RNA Library Prep and Target Enrichment workflow. See the [Sample Preparation](#) chapter for complete instructions for use of the VWorks protocols for sample processing.

Table 7 Overview of VWorks protocols and runsets used during the workflow

Workflow Step	VWorks Protocols Used for Agilent NGS Workstation automation
<ul style="list-style-type: none"> Purify poly(A) RNA using oligo(dT) beads Chemically fragment the poly(A) RNA Synthesize first-strand cDNA 	mRNA_Purification_v1.0.pro
Purify first-strand cDNA using AMPure XP beads	AMPureXP_v1.1.pro:First Strand
<ul style="list-style-type: none"> Synthesize second-strand cDNA Repair DNA ends Purify end-repaired DNA dA-tail DNA 3'-ends Ligate adaptors Purify adaptor-ligated DNA 	LibraryPrep_RNASeq_ILM_v1.1.rst
Amplify adaptor-ligated cDNA library	Pre-CapturePCR_RNASeq_ILM_v1.0.pro
Purify library amplicons using AMPure XP beads	AMPureXP_v1.1.pro:Pre-Capture PCR
Aliquot 100 ng of prepped libraries for hybridization	Aliquot_Libraries_v1.0.pro
Hybridize prepped DNA to Capture Library	SureSelectHybridization_v1.0.pro
Capture and wash DNA hybrids	SureSelectCapture&Wash_v1.0.rst
Add index tags by PCR	Post-CapturePCR_RNASeq_ILM_v1.0.pro
Purify indexed amplicons using AMPure XP beads	AMPureXP_v1.1.pro:Post-Capture PCR

Experimental Setup Considerations for Automated Runs

Agilent SureSelect Automated Strand-Specific RNA Library Prep runs may include 1, 2, 3, 4, 6, or 12 columns (equivalent to 8, 16, 24, 32, 48, or 96 wells) of RNA samples to be prepared for sequencing on the Illumina platform. Plan your experiments using complete columns of samples.

Table 8 Columns to Samples Equivalency

Number of Columns Processed	Total Number of Samples Processed
1	8
2	16
3	24
4	32
6	48
12	96

The number of columns or samples that may be processed using the supplied reagents will depend on the experimental design. For greatest efficiency of reagent use, plan experiments using at least 3 columns per run. Each 96-reaction kit contains sufficient reagents for 96 reactions configured as 4 runs of 3 columns of samples per run.

Considerations for Placement of RNA Samples in 96-well Plates for Automated Processing

- The Agilent NGS Workstation processes samples column-wise beginning at column 1. RNA samples should be loaded into 96-well plates column-wise, in well order A1 to H1, then A2 to H2, ending with A12 to H12. When processing partial runs with <12 sample columns, do not leave empty columns between sample columns; always load the plate using the left-most column that is available.
- For sample indexing by PCR (see [Figure 2](#)), you will need to prepare a separate plate containing the indexing primers. Assign the wells to be indexed with their respective indexing primers during experimental design.

Considerations for Equipment Setup

- Some workflow steps require the rapid transfer of sample plates between the Bravo deck and a thermal cycler. Locate your thermal cycler in close proximity to the Agilent NGS Workstation to allow rapid and efficient plate transfer.
- Several workflow steps require that the sample plate be sealed using the PlateLoc thermal microplate sealer on the Agilent NGS Workstation, and then centrifuged to collect any dispersed liquid. To maximize efficiency, locate the centrifuge in close proximity to the Agilent NGS Workstation.

PCR Plate Type Considerations

Automation protocols include several liquid-handling steps in which reagents are dispensed to PCR plates in preparation for transfer to a thermal cycler. For these steps you must specify the PCR plate type to be used on the SureSelect_RNA_ILM.VWForm to allow correct configuration of the liquid handling components for the PCR plate type. Before you begin the automation protocol, make sure that you are using a supported PCR plate type. The PCR plate type to be used in the protocol is specified using the menu below. Vendor and part number information is provided for the supported plate types in [Table 9](#).

2) Select PCR Plate labware for Thermal Cycling

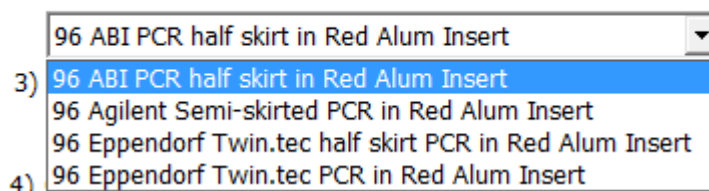


Table 9 Ordering information for supported PCR plates

Description in VWorks menu	Vendor and part number
96 ABI PCR half-skirted plates (MicroAmp Optical plates)	Thermo Fisher Scientific p/n N8010560
96 Agilent semi-skirted PCR plate	Agilent p/n 401334
96 Eppendorf Twin.tec half-skirted PCR plates	Eppendorf p/n 951020303
96 Eppendorf Twin.tec PCR plates (full-skirted)	Eppendorf p/n 951020401 or 951020619

CAUTION

The plates listed in [Table 9](#) are compatible with the Agilent NGS Bravo and associated VWorks automation protocols, designed to support use of various thermal cyclers.

Accordingly, some plates listed in [Table 9](#) are not compatible with the recommended SureCycler 8800 Thermal Cycler. When using the SureCycler 8800 Thermal Cycler in the SureSelect automation workflow, use 96 Agilent semi-skirted PCR plates.

When using a different thermal cycler in the workflow, be sure to select a PCR plate that is compatible with your thermal cycler and that is listed in [Table 9](#).



3 Sample Preparation

- Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA 34
- Step 2. Purify first-strand cDNA using AMPure XP beads 55
- Step 3. Prepare cDNA libraries for Hybridization 59
- Step 4. Amplify cDNA libraries by PCR 67
- Step 5. Purify amplified DNA using AMPure XP beads 75
- Step 6. Assess library DNA quantity and quality 79

This section contains instructions for RNA library preparation specific to the Illumina paired-read sequencing platform and to automated processing using the Agilent NGS Workstation.

For each sample to be sequenced, individual library preparations are performed in separate wells of a 96-well plate. The samples are then target-enriched and indexed by PCR amplification allowing multiplexing of up to 96 samples for sequencing on Illumina platforms.



3 Sample Preparation

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

In this step, automation protocol mRNA_Purification_v1.0.pro is used to complete multiple steps of the RNA Library Preparation workflow. First, poly(A) RNA is purified from total RNA using two serial rounds of binding to oligo(dT) magnetic particles. After purification, the poly(A) RNA is chemically-fragmented to the appropriate size and then is converted to first-strand cDNA.

Total RNA samples containing 200 ng to 4 µg RNA are suitable for the mRNA library preparation automation protocol. Each total RNA sample must be prepared for the run in 25 µL of nuclease-free water.

NOTE

For optimal performance, total RNA samples should have an RNA Integrity Number (RIN) of 8 or more, based on analysis using Agilent's 2100 Bioanalyzer.

Consider preparing an additional sequencing library in parallel, using a high-quality control RNA sample, such as Agilent QPCR Human Reference Total RNA (p/n 750500). Use of this control is especially recommended during the first run of the protocol, to verify that all protocol steps are being successfully performed. Routine use of this control is helpful for any required troubleshooting, in order to differentiate any performance issues related to RNA input from other factors.

A workstation operator must be present during this automation protocol to transfer plates between the workstation, which completes most liquid handling steps, and the thermal cycler, which is used for several incubation steps. In addition, the operator must prepare and dispense a master mix immediately before it is used in the automation protocol (see [step 30](#) on [page 50](#)).

Prepare the workstation

- 1 Open the SureSelect setup form using the SureSelect_RNA_ILM.VWForm shortcut on your desktop.
- 2 Log in to the VWorks software.

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

- 3 Turn on the ThermoCube, set to 0°C, at position 9 of the Bravo deck. Be sure that the chiller reservoir contains at least 300 mL of 25% ethanol.
- 4 Clear the Labware MiniHub and BenchCel of all plates and tip boxes.

Prepare reagents for the run

- 5 Bring the reagents listed in [Table 10](#) to room temperature and thaw the reagents listed in [Table 11](#) on ice.

Table 10 Reagents brought to room temperature before use in protocol

Storage Location	Kit Component	Where Used in Protocol
SureSelect Poly-A Selection Module [*] , 4°C	Oligo(dT) Microparticles (bottle)	page 37
	Bead Washing Buffer [†] (bottle)	page 37
	Bead Elution Buffer [‡] (bottle)	page 37
	Bead Binding Buffer ^{**} (bottle)	page 37

* May also be labeled as *SureSelect Strand Specific RNA Library Prep, ILM, Box 2*.

† May also be labeled as *RNA Seq Bead Washing Buffer*.

‡ May also be labeled as *RNA Seq Bead Elution Buffer*.

** May also be labeled as *RNA Seq Bead Binding Buffer*.

Table 11 Reagents thawed and held on ice before use in protocol

Storage Location	Kit Component	Where Used in Protocol
SureSelect RNA Library Prep, ILM (Pre PCR) [†] , -20°C	Fragmentation Mix [†] (bottle)	page 38
	First Strand Master Mix [‡] (tube with orange cap)	page 51

* May also be labeled as *SureSelect Strand Specific RNA Library Prep, ILM, Box 1*.

† May also be labeled as *RNA Seq Fragmentation Mix*.

‡ May also be labeled as *RNA Seq First Strand Master Mix*.

3 Sample Preparation

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

- 6 Locate or prepare a stock solution of 4 µg/µL Actinomycin D in DMSO. When first prepared, aliquot the stock solution into single-use volumes (typically 3 µL) and store the aliquots at -20°C, protected from light. Do not subject the aliquots to multiple freeze-thaw cycles. The aliquots may be stored for up to one year before use in the library preparation protocol.

A 3-µL aliquot of this DMSO stock solution will be used on [page 50](#) to prepare a fresh dilution of 120 ng/µL Actinomycin D in water for the run.

CAUTION

To ensure strand specificity, take care to follow the storage instructions above for the 4 µg/µL Actinomycin D in DMSO stock solution, including the one-year maximum storage duration.

Prepare the RNA samples source plate

- 7 Place 25 µL of each RNA sample (0.2–4 µg RNA in nuclease-free water) into the wells of a 96-well Eppendorf twin.tec plate. Load samples into the plate column-wise, in well order A1 to H1, then A2 to H2, ending with A12 to H12 for processing on the Agilent NGS Workstation.

NOTE

SureSelect Strand-Specific RNA Library Prep runs may include 1, 2, 3, 4, 6, or 12 columns of the plate. See [Using the Agilent NGS Workstation for SureSelect RNA Library Preparation](#) for additional sample placement considerations.

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

Prepare the oligo(dT) beads and mRNA purification source plates

When preparing each of the source plates below, add the indicated amount of reagent to wells of the source plate corresponding to the total RNA sample wells in [step 7](#) above. For example, for 3-column runs, fill source well plate wells A1 to H3, but leave wells A4 to H12 empty.

- 8** Prepare the oligo(dT) beads source plate.
 - a** Vortex the Oligo(dT) Microparticles until the suspension appears homogeneous and consistent in color. If bead aggregates are still present after vortexing, mix thoroughly by pipetting up and down until the suspension appears homogeneous.
 - b** In a PCR plate that is compatible with the thermal cycler to be used in the run, place 25 μL of the homogeneous Oligo(dT) bead suspension into each well to be used for sample purification.
- 9** Prepare the Bead Binding Buffer source plate. Place 30 μL of Bead Binding Buffer into wells of a 96-well Eppendorf twin.tec plate. Fill each well that corresponds to an RNA sample well.
- 10** Prepare the Bead Elution Buffer source plate. Place 30 μL of Bead Elution Buffer into wells of a 96-well Eppendorf twin.tec plate. Fill each well that corresponds to an RNA sample well.
- 11** Prepare the Bead Wash Buffer source plate. Place 410 μL of Bead Washing Buffer into wells of a Nunc DeepWell plate. Fill each well that corresponds to an RNA sample well.

3 Sample Preparation

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

Prepare the master mix source plate

12 Prepare the master mix source plate by adding the appropriate volume of Fragmentation Mix (see Table 12) to all wells of Column 1 of a Nunc DeepWell plate. The configuration of the source plate is shown in Figure 3.

Table 12 Preparation of the Master Mix Source Plate for mRNA_Purification_v1.0.pro

Master Mix Solution	Position on Source Plate	Volume of Master Mix added per Well of Nunc Deep Well Source Plate					
		1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
Fragmentation Mix	Column 1 (A1-H1)	28.5 µL	47.5 µL	66.5 µL	85.5 µL	123.5 µL	247.0 µL

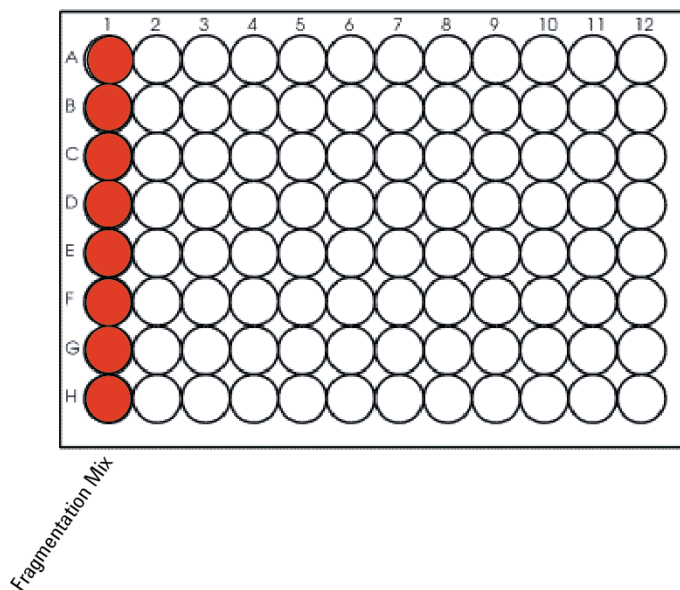


Figure 3 Initial configuration of master mix source plate for mRNA_Purification_v1.0.pro

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

Load the Agilent NGS Workstation

13 Load the Labware MiniHub according to [Table 13](#), using the plate orientations shown in [Figure 4](#).

Table 13 Initial MiniHub configuration for mRNA_Purification_v1.0.pro

Vertical Shelf Position	Cassette 1	Cassette 2	Cassette 3	Cassette 4
Shelf 5 (Top)	Empty Nunc DeepWell plate	Empty Nunc DeepWell plate	Empty Nunc DeepWell plate	Empty
Shelf 4	Empty	Bead Binding Buffer in twin.tec plate	Empty	Empty
Shelf 3	Empty	Bead Elution Buffer in twin.tec plate	Empty	Empty
Shelf 2	Empty tip box	Empty	Bead Wash Buffer in Nunc DeepWell plate	Empty
Shelf 1 (Bottom)	New tip box	Empty	Empty	Empty tip box

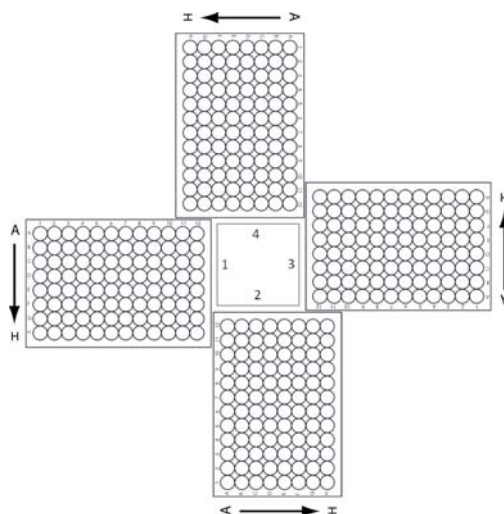


Figure 4 Agilent Labware MiniHub plate orientation. For Thermo Scientific reservoirs, place the notched corner facing the center of the hub.

3 Sample Preparation

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

14 Load the Bravo deck according to [Table 14](#).

Table 14 Initial Bravo deck configuration for mRNA_Purification_v1.0.pro

Location	Content
1	Empty waste reservoir (Axygen 96 Deep Well Plate, square wells)
4	Oligo(dT) beads in PCR plate seated on red insert (PCR plate type must be specified on setup form under step 2)
6	Empty PCR plate seated on red insert (PCR plate type must be specified on setup form under step 2)
7	Total RNA samples in twin.tec plate
9	Master Mix Source Plate seated on silver insert (Nunc DeepWell; see Figure 3 on page 38 for column content)

15 Load the BenchCel Microplate Handling Workstation according to [Table 15](#).

Table 15 Initial BenchCel configuration for mRNA_Purification_v1.0.pro

No. of Columns Processed	Rack 1	Rack 2	Rack 3	Rack 4
1	1 Tip box	Empty	Empty	Empty
2	2 Tip boxes	Empty	Empty	Empty
3	3 Tip boxes	Empty	Empty	Empty
4	3 Tip boxes	Empty	Empty	Empty
6	5 Tip boxes	Empty	Empty	Empty
12	9 Tip boxes	Empty	Empty	Empty

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

Run VWorks protocol mRNA_Purification_v1.0.pro

16 On the SureSelect setup form, under **Select Protocol to Run**, select **mRNA_Purification_v1.0.pro**.

17 Under **Select PCR plate labware for Thermal Cycling**, select the specific type of PCR plate that was loaded on Bravo deck positions 4 and 6. The plate type selected must be compatible with the thermal cycler to be used for incubation steps during the protocol.

NOTE

During run setup, be sure to use the plate type selected from this menu at positions 4 and 6 of the Bravo deck. In addition, when the workstation issues prompts to add plates to position 4 or 6 during the run, use only the same PCR plate type specified here.

18 Select the number of columns of samples to be processed. Runs must include 1, 2, 3, 4, 6, or 12 columns.

3) Select Number of Columns of Samples

4) 1
2
3
4
6
12

Display Initial Workstation Setup

Clear Setup

19 Click **Display Initial Workstation Setup**.



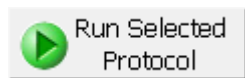
20 Verify that the NGS workstation has been set up as displayed in the Workstation Setup region of the form.

	MiniHub Cassette 1	MiniHub Cassette 2	MiniHub Cassette 3	MiniHub Cassette 4
Shelf 5				
Shelf 4				

3 Sample Preparation

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

21 When verification is complete, click **Run Selected Protocol**.



NOTE

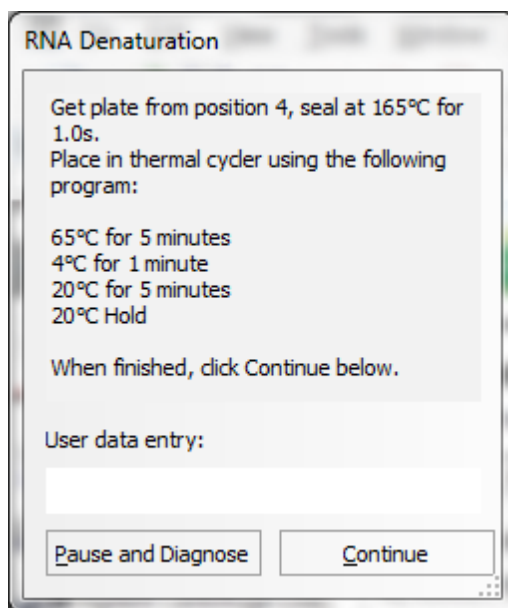
If workstation devices do not respond when you start the run, but activity is recorded in the Log, verify that VWorks is not running in Simulation mode. See [page 27](#) for more information.

Running the mRNA_Purification_v1.0.pro protocol takes approximately 90 minutes, including four incubation periods on the thermal cycler.

During the automation protocol run, a workstation operator must be present to transfer plates between the workstation and thermal cycler when prompted, as detailed on the following pages.

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

22 When the workstation has finished combining the RNA samples with the oligo(dT) beads, you will be prompted by VWorks as shown below.



- a Remove the plate from position 4 of the Bravo deck and seal the plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 1.0 seconds.
- b Briefly spin the plate in a centrifuge or mini-plate spinner to collect the liquid without pelleting the beads.
- c Transfer the PCR plate to a thermal cycler (with the heated lid ON) and run the RNA denaturation + bead binding program shown in [Table 16](#). After transferring the plate, click **Continue** on the VWorks screen.

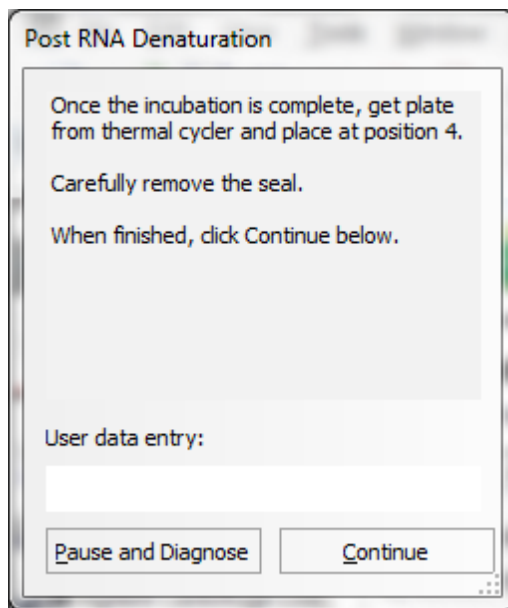
Table 16 Thermal cycler program for RNA denaturation and RNA-bead binding

Step	Temperature	Time
Step 1	65°C	5 minutes
Step 2	4°C	1 minute
Step 3	20°C	5 minutes
Step 4	20°C	Hold

3 Sample Preparation

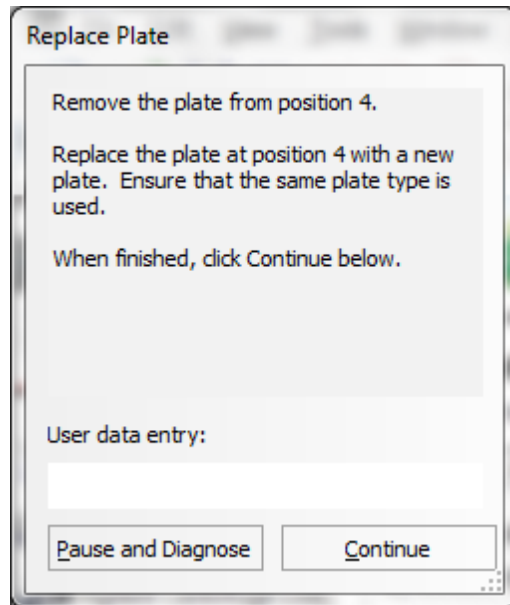
Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

- 23** After the thermal cycler reaches the 20°C Hold step, and when prompted by the dialog below, transfer the RNA sample plate to position 4 of the Bravo deck, seated in the red insert. Carefully unseal the plate, then click **Continue**.



Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

- 24 When the workstation has finished collecting and washing the bead-bound RNA samples, you will be prompted by VWorks as shown below.

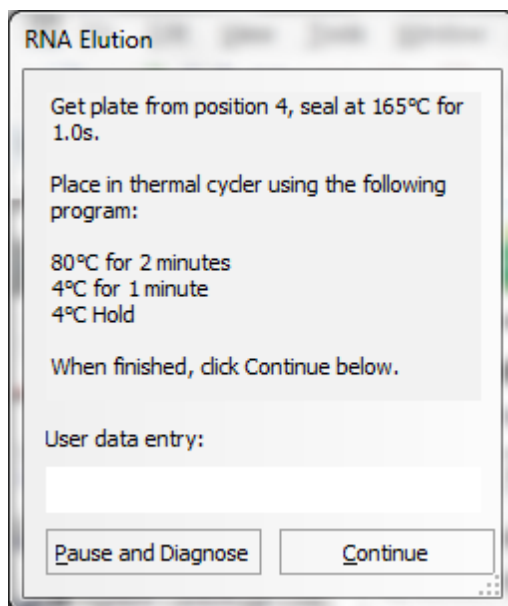


- a Remove and discard the PCR plate from position 4 of the Bravo deck.
- b Place a fresh PCR plate at position 4, seated in the red insert. The PCR plate type added here must be the same plate type as the one removed and as was specified during the run setup.
- c After positioning the plate, click **Continue** on the VWorks screen.

3 Sample Preparation

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

25 When the workstation has finished collecting and washing the bead-bound RNA samples, you will be prompted to transfer the plate to the thermal cycler for the RNA Elution step as shown below.



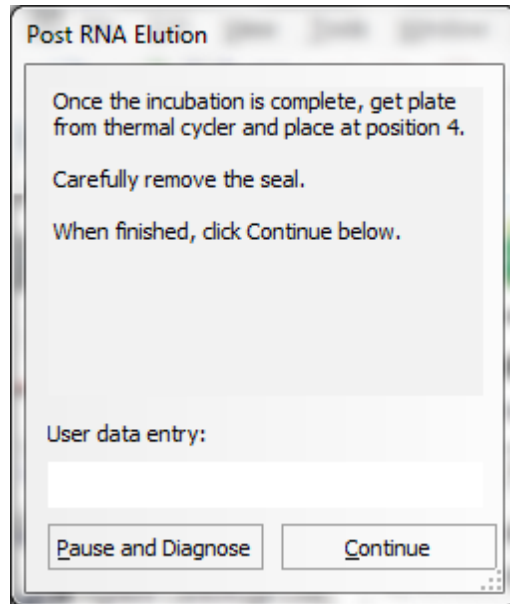
- a** Remove the plate from position 4 of the Bravo deck and seal the plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 1.0 seconds.
- b** Briefly spin the plate in a centrifuge or mini-plate spinner to collect the liquid without pelleting the beads.
- c** Transfer the PCR plate to a thermal cycler (with the heated lid ON) and run the RNA elution program shown in [Table 17](#). After transferring the plate, click **Continue** on the VWorks screen.

Table 17 Thermal cycler program for RNA elution

Step	Temperature	Time
Step 1	80°C	2 minutes
Step 2	4°C	1 minute
Step 3	4°C	Hold

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

- 26** After the thermal cycler reaches the 4°C Hold step and when prompted by the dialog below, transfer the RNA sample plate to position 4 of the Bravo deck, seated in the red insert. Carefully unseal the plate, then click **Continue**.

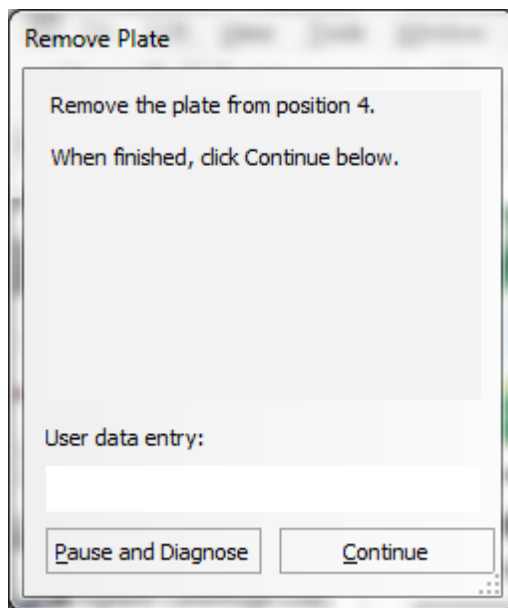


- 27** The workstation adds Bead Binding Buffer to the eluted RNA samples and then holds the samples at room temperature for 5 minutes to allow the poly(A) RNA to re-bind the beads.

3 Sample Preparation

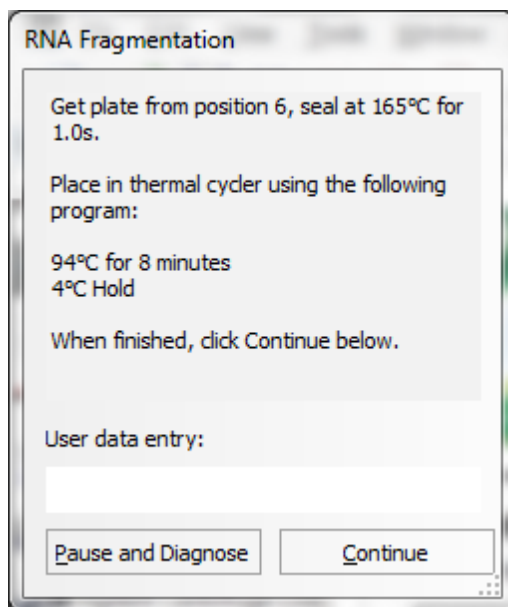
Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

- 28** When the workstation has finished the collecting and washing the bound RNA samples in this second round of purification, you will be prompted by VWorks as shown below. Remove and discard the PCR plate from position 4, then click **Continue**.



Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

- 29** The workstation adds Fragmentation Mix to the bead-bound RNA samples in preparation for the RNA fragmentation step. When the workstation has finished, you will be prompted by VWorks as shown below.



- a** Remove the plate from position 6 of the Bravo deck and seal the plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 1.0 seconds.
- b** Briefly spin the plate in a centrifuge or mini-plate spinner to collect the liquid.
- c** Transfer the PCR plate to a thermal cycler (with the heated lid ON) and run the RNA fragmentation program shown in [Table 18](#). After transferring the plate, click **Continue** on the VWorks screen.

Table 18 Thermal cycler program for RNA fragmentation

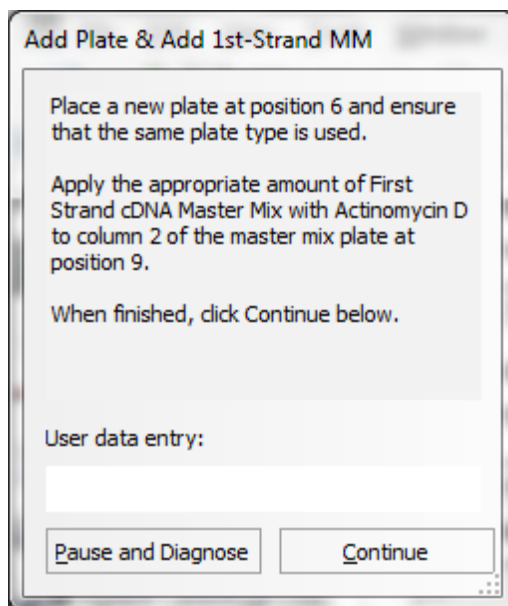
Step	Temperature	Time
Step 1	94°C	8 minutes
Step 2	4°C	Hold

3 Sample Preparation

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

30 During the 8-minute incubation step, prepare the reagents and workstation for first-strand cDNA synthesis:

- a** When prompted by the dialog below, place a fresh PCR plate (use the plate type specified during the run setup) at position 6, seated in the red insert. Proceed immediately to [step b](#), below.



- b** Prepare a fresh 120 ng/ μ L Actinomycin D dilution in water from a stock solution of 4 μ g/ μ L Actinomycin D in DMSO, according to [Table 19](#).

CAUTION

To ensure strand-specificity, you must prepare the 120 ng/ μ L Actinomycin D dilution in water immediately before use in [step d](#), below.

Table 19 Preparation of 120 ng/ μ L Actinomycin D

Reagent	Volume for up to 12-column run (includes excess)
Actinomycin D (4 μ g/ μ L in DMSO)	3 μ L
Nuclease-free water	97 μ L
Total	100 μL

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

- c Vortex the thawed vial of First Strand Master Mix 5 seconds at high speed to ensure homogeneity.

CAUTION

SureSelect RNA Library Prep master mixes are viscous. Mix thoroughly by vortexing at high speed for 5 seconds before removing an aliquot for use and after combining the master mix with other solutions. Pipetting up and down is not sufficient to mix this reagent.

- d Prepare the appropriate amount of First Strand Master Mix + Actinomycin D mixture, on ice, according to the table below. Mix by vortexing at high speed for 5 seconds, then spin briefly and keep on ice.

Table 20 Preparation of First Strand Master Mix/Actinomycin D mixture

Reagent	Volume for 1 Library	Volume for 1 Column	Volume for 2 Columns	Volume for 3 Columns	Volume for 4 Columns	Volume for 6 Columns	Volume for 12 Columns
First Strand Master Mix	8.0 µL	98.4 µL	196.8 µL	262.4 µL	360.8 µL	492.0 µL	918.4 µL
Actinomycin D (120 ng/µL in H ₂ O)	0.5 µL	6.2 µL	12.3 µL	16.4 µL	22.6 µL	30.8 µL	57.4 µL
Total Volume	8.5 µL	104.6 µL	209.1 µL	278.8 µL	383.4 µL	522.8 µL	975.8 µL

- e Add the volume listed in [Table 21](#) of the First Strand Master Mix + Actinomycin D mixture to column 2 of the Master Mix source plate at position 9 of the Bravo deck. The final configuration of the source plate is shown in [Figure 6](#). After adding the master mix to the source plate, click **Continue** on the VWorks screen.

Table 21 Preparation of the Master Mix Source Plate for mRNA_Purification_v1.0.pro

Master Mix Solution	Position on Source Plate	Volume of Master Mix added per Well of Nunc Deep Well Source Plate					
		1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
First Strand Master Mix + Actinomycin D mixture	Column 2 (A2-H2)	12.0 µL	25.1 µL	33.8 µL	46.9 µL	64.3 µL	120.9 µL

3 Sample Preparation

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

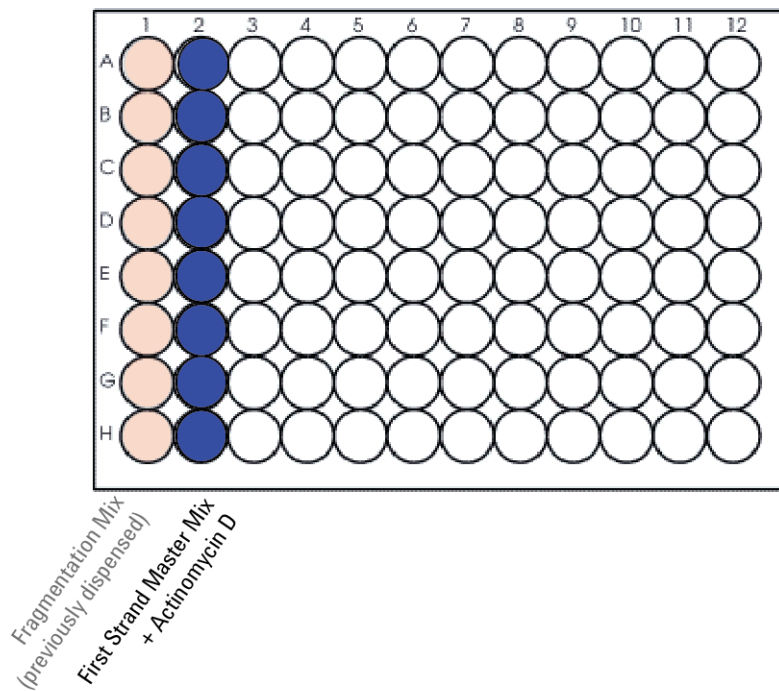
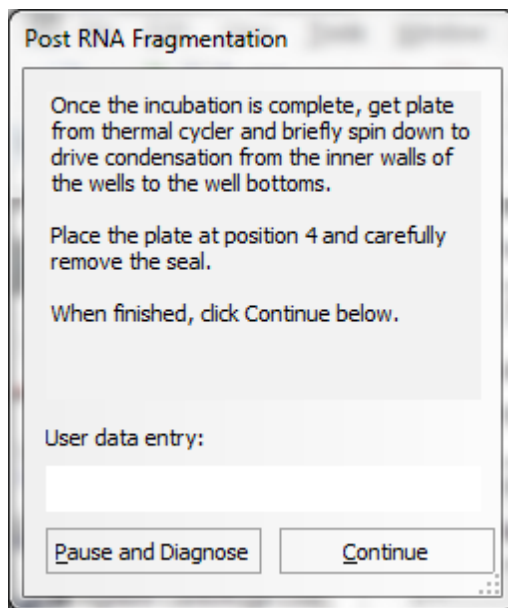


Figure 5 Final configuration of master mix source plate for mRNA_Purification_v1.0.pro

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

- 31** With the RNA sample plate still on the thermal cycler, the workstation prepares the remaining components for first-strand cDNA synthesis. When the workstation has finished, you will be prompted by VWorks as shown below.

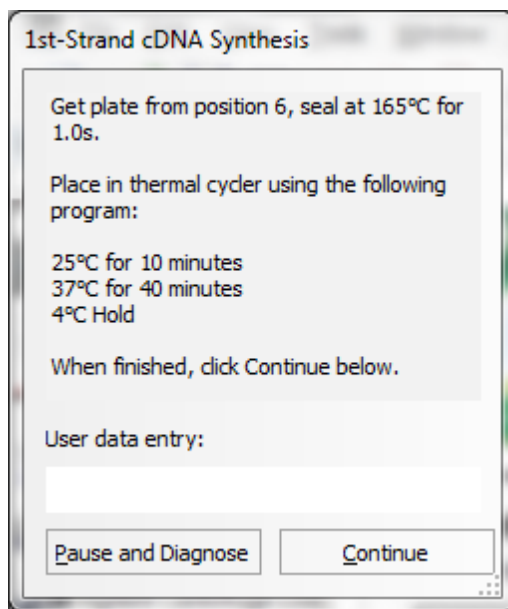


- a** After the thermal cycler reaches the 4°C Hold step for the RNA fragmentation program (Table 18), remove the plate from the thermal cycler and briefly spin in a centrifuge or mini-plate spinner to collect the liquid.
- b** Place the RNA sample plate on position 4 of the Bravo deck, seated in the red insert.
- c** Carefully unseal the plate, then click **Continue**.

3 Sample Preparation

Step 1. Purify poly(A) RNA/fragment RNA/synthesize first strand cDNA

- 32** The workstation removes the fragmented RNA samples from the bead-containing wells and combines the samples with First Strand Master Mix + Actinomycin D. When the workstation has finished, you will be prompted by VWorks as shown below.



- a** Remove the plate from position 6 of the Bravo deck and seal the plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 1.0 seconds.
- b** Briefly spin the plate in a centrifuge or mini-plate spinner to collect the liquid.
- c** Transfer the PCR plate to a thermal cycler (with the heated lid ON) and run the first-strand cDNA synthesis program shown in [Table 22](#). After transferring the plate, click **Continue** on the VWorks screen.

Table 22 Thermal cycler program for first-strand cDNA synthesis

Step	Temperature	Time
Step 1	25°C	10 minutes
Step 2	37°C	40 minutes
Step 3	4°C	Hold

Step 2. Purify first-strand cDNA using AMPure XP beads

In this step, the Agilent NGS Workstation transfers AMPure XP beads and first-strand cDNA samples to a Nunc DeepWell plate and then collects and washes the bead-bound DNA.

Prepare the workstation and reagents

- 1 Leave tip boxes on shelves 1 and 2 in cassette 1 of the Labware MiniHub from the previous mRNA_Purification_v1.0.pro run. Otherwise, clear the remaining positions of the MiniHub and BenchCel of plates and tip boxes.
- 2 Gently wipe down the Labware MiniHub, Bravo deck, and BenchCel with a DNA Away decontamination wipe.
- 3 Turn on the ThermoCube, set to 0°C, at position 9 of the Bravo deck. Be sure that the chiller reservoir contains at least 300 mL of 25% ethanol.
- 4 Let the AMPure XP beads come to room temperature for at least 30 minutes. *Do not freeze the beads at any time.*
- 5 Mix the bead suspension well so that the reagent appears homogeneous and consistent in color.
- 6 Prepare a Nunc DeepWell source plate for the beads by adding 51 µL of homogeneous AMPure XP beads per well, for each well to be processed.
- 7 Prepare a Thermo Scientific reservoir containing 15 mL of nuclease-free water.
- 8 Prepare a separate Thermo Scientific reservoir containing 45 mL of freshly-prepared 70% ethanol.

3 Sample Preparation

Step 2. Purify first-strand cDNA using AMPure XP beads

- 9 Load the Labware MiniHub according to [Table 23](#), using the plate orientations shown in [Figure 4](#).

Table 23 Initial MiniHub configuration for AMPureXP_v1.1.pro:First Strand

Vertical Shelf Position	Cassette 1	Cassette 2	Cassette 3	Cassette 4
Shelf 5 (Top)	Empty Nunc DeepWell plate	Empty	Empty	Empty
Shelf 4	Empty	Empty	Empty	Empty
Shelf 3	Empty	Empty Eppendorf twin.tec Plate	Empty	Empty
Shelf 2	Waste tip box (retained from mRNA_Purification protocol)*	Nuclease-free water reservoir from step 7	AMPure XP beads in Nunc DeepWell plate from step 6	Empty
Shelf 1 (Bottom)	Clean tip box (retained from mRNA_Purification protocol)*	70% ethanol reservoir from step 8	Empty	Empty tip box

* The tip boxes retained in Cassette 1 are not shown on the VWorks Workstation Setup table. These tip boxes are not used in AMPureXP_v1.1.pro:First Strand but are used in a later protocol. This labware should be retained in the MiniHub to ensure that empty and full tip positions are properly defined for the subsequent protocol.

- 10 Load the Bravo deck according to [Table 24](#).

Table 24 Initial Bravo deck configuration for AMPureXP_v1.1.pro:First Strand

Location	Content
1	Empty waste reservoir (Axygen 96 Deep Well Plate, square wells)
9	First-strand cDNA samples in PCR plate seated on red insert (PCR plate type must be specified on setup form under step 2)

Step 2. Purify first-strand cDNA using AMPure XP beads

11 Load the BenchCel Microplate Handling Workstation according to Table 25.

Table 25 Initial BenchCel configuration for AMPureXP_v1.1.pro:First Strand

No. of Columns Processed	Rack 1	Rack 2	Rack 3	Rack 4
1	1 Tip box	Empty	Empty	Empty
2	1 Tip box	Empty	Empty	Empty
3	2 Tip boxes	Empty	Empty	Empty
4	2 Tip boxes	Empty	Empty	Empty
6	3 Tip boxes	Empty	Empty	Empty
12	6 Tip boxes	Empty	Empty	Empty

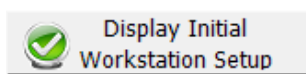
Run VWorks protocol AMPureXP_v1.1.pro:First Strand

12 On the SureSelect setup form, under **Select Protocol to Run**, select **AMPureXP_v1.1.pro:First Strand**.

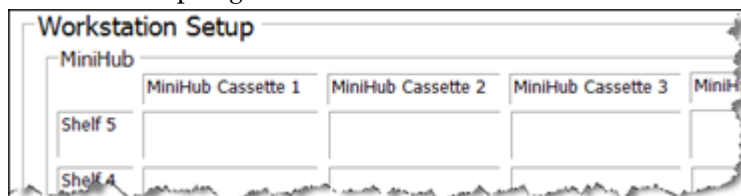
13 Under **Select PCR plate labware for Thermal Cycling**, select the specific type of PCR plate containing the cDNA samples at position 9.

14 Select the number of columns of samples to be processed. Runs must include 1, 2, 3, 4, 6, or 12 columns.

15 Click **Display Initial Workstation Setup**.



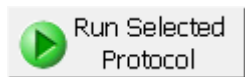
16 Verify that the NGS workstation has been set up as displayed in the Workstation Setup region of the form.



3 Sample Preparation

Step 2. Purify first-strand cDNA using AMPure XP beads

17 When verification is complete, click **Run Selected Protocol**.



Running the AMPureXP_v1.1.pro:First Strand protocol takes approximately 45 minutes. During this time, you can prepare the purification reagents for the Library Prep automation protocol, as described on [page 60](#).

Once the AMPureXP_v1.1.pro:First Strand protocol is complete, the purified cDNA samples are located in the Eppendorf plate at position 7 of the Bravo deck. Proceed immediately to “[Step 3. Prepare cDNA libraries for Hybridization](#)” on page 59.

Step 3. Prepare cDNA libraries for Hybridization

This step is automated using the LibraryPrep_RNASeq_ILM_v1.1.rst runset. During the runset, the Agilent NGS Workstation completes second-strand cDNA library synthesis and end modification steps, including end-repair, A-tailing, and adaptor ligation. After certain modification steps, the Agilent NGS Workstation purifies the prepared cDNA using AMPure XP beads.

This step uses the SureSelect^{XT} RNA Reagent Kit components listed in Table 26 in addition to the purification reagents prepared for use on page 60. Thaw each reagent vial and keep on ice. Vortex each vial for 5 seconds at high speed to mix before use.

Table 26 Reagents for automation runset LibraryPrep_RNASeq_ILM_v1.1.rst

Storage Location	Kit Component	Where Used in Protocol
SureSelect RNA Library Prep, ILM (Pre PCR) [*] , -20°C	Second Strand Enzyme Mix [†] (bottle)	page 61
	Second Strand Oligo Mix [‡] (tube with yellow cap)	page 61
	SureSelect Ligation Master Mix (tube with purple cap)	page 61
	SureSelect Oligo Adaptor Mix (tube with blue cap)	page 61
	dA Tailing Master Mix ^{**} (bottle)	page 62

* May also be labeled as *SureSelect Strand Specific RNA Library Prep, ILM, Box 1*.

† May also be labeled as *RNA Seq Second Strand + End Repair Enzyme Mix*.

‡ May also be labeled as *RNA Seq Second Strand + End Repair Oligo Mix*.

** May also be labeled as *RNA Seq dA Tailing Master Mix*.

Prepare the workstation

- 1 Turn on the ThermoCube, set to 0°C, at position 9 of the Bravo deck. Be sure that the chiller reservoir contains at least 300 mL of 25% ethanol.
- 2 Leave tip boxes on shelves 1 and 2 in cassette 1 of the Labware MiniHub from the previous mRNA_Purification_v1.0.pro run. Otherwise, clear the remaining MiniHub and BenchCel positions of plates and tip boxes.

3 Sample Preparation

Step 3. Prepare cDNA libraries for Hybridization

- 3 Pre-set the temperature of Bravo deck position 4 to 14°C and of position 6 to 4°C using the Inheco Multi TEC control touchscreen, as described in [Setting the Temperature of Bravo Deck Heat Blocks](#). On the control touchscreen, Bravo deck positions 4 corresponds to CPAC 2, position 1, while deck position 6 corresponds to CPAC 2, position 2.

Prepare the purification reagents

- 4 Verify that the AMPure XP bead suspension is at room temperature. *Do not freeze the beads at any time.*
- 5 Mix the bead suspension well so that the reagent appears homogeneous and consistent in color.
- 6 Prepare a Nunc DeepWell source plate for the beads by adding 160 µL of homogeneous AMPure XP beads per well, for each well to be processed.
- 7 Prepare a Thermo Scientific reservoir containing 20 mL of nuclease-free water.
- 8 Prepare a separate Thermo Scientific reservoir containing 150 mL of freshly-prepared 70% ethanol.

Prepare the master mix source plate

- 9 Vortex the thawed vials of Second Strand Enzyme Mix, SureSelect Ligation Master Mix, and dA Tailing Master Mix for 5 seconds at high speed to ensure homogeneity.

CAUTION

SureSelect RNA Library Prep master mixes are viscous. Mix thoroughly by vortexing at high speed for 5 seconds before removing an aliquot for use and after combining the master mix with other solutions. Pipetting up and down is not sufficient to mix these reagents.

- 10 Prepare the appropriate amount of Second Strand + End Repair Master Mix according to [Table 27](#) below.

Table 27 Preparation of Second Strand + End Repair Master Mix for LibraryPrep_RNASeq_ILM_v1.1.rst

Reagent	Volume for 1 Library	Volume for 1 Column	Volume for 2 Columns	Volume for 3 Columns	Volume for 4 Columns	Volume for 6 Columns	Volume for 12 Columns
Second Strand Enzyme Mix	25.0 µL	307.5 µL	615 µL	820 µL	1127.5 µL	1640 µL	3075 µL
Second Strand Oligo Mix	5.0 µL	61.5 µL	123 µL	164 µL	225.5 µL	328 µL	615 µL
Total Volume	30 µL	369 µL	738 µL	984 µL	1353 µL	1968 µL	3690 µL

- 11 Prepare the appropriate amount of Adaptor Ligation Master Mix, containing the SureSelect Ligation Master Mix and the adaptors, according to [Table 28](#) below.

Table 28 Preparation of Adaptor Ligation Master Mix for LibraryPrep_RNASeq_ILM_v1.1.rst

Reagent	Volume for 1 Library	Volume for 1 Column	Volume for 2 Columns	Volume for 3 Columns	Volume for 4 Columns	Volume for 6 Columns	Volume for 12 Columns
Nuclease-free water	2.5 µL	30.8 µL	61.5 µL	82.0 µL	112.8 µL	164.0 µL	307.5 µL
SureSelect Ligation Master Mix	5.0 µL	61.5 µL	123.0 µL	164.0 µL	225.5 µL	328.0 µL	615.0 µL
SureSelect Oligo Adaptor Mix	5.0 µL	61.5 µL	123.0 µL	164.0 µL	225.5 µL	328.0 µL	615.0 µL
Total Volume	12.5 µL	153.8 µL	307.5 µL	410.0 µL	563.8 µL	820.0 µL	1537.5 µL

3 Sample Preparation

Step 3. Prepare cDNA libraries for Hybridization

12 Using the same Nunc DeepWell master mix source plate that was used for the mRNA_Purification_v1.0.pro run, prepare the Library Prep master mix source plate. Add the volumes indicated in [Table 29](#) of each master mix to all wells of the indicated column of the plate. Keep the master mixes on ice during the aliquoting steps. The final configuration of the master mix source plate is shown in [Figure 6](#).

Table 29 Preparation of the Master Mix Source Plate for LibraryPrep_RNASeq_ILM_v1.1.rst

Master Mix Solution	Position on Source Plate	Volume of Master Mix added per Well of Nunc Deep Well Source Plate					
		1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
Second Strand + End Repair Master Mix (from Table 27)	Column 3 (A3-H3)	42.4 μ L	88.5 μ L	119.3 μ L	165.4 μ L	242.3 μ L	457.5 μ L
dA Tailing Master Mix	Column 4 (A4-H4)	30.0 μ L	50.0 μ L	70.0 μ L	90.0 μ L	130.0 μ L	260.0 μ L
Adaptor Ligation Master Mix (from Table 28)	Column 5 (A5-H5)	17.7 μ L	36.9 μ L	49.7 μ L	68.9 μ L	100.9 μ L	190.6 μ L

Step 3. Prepare cDNA libraries for Hybridization

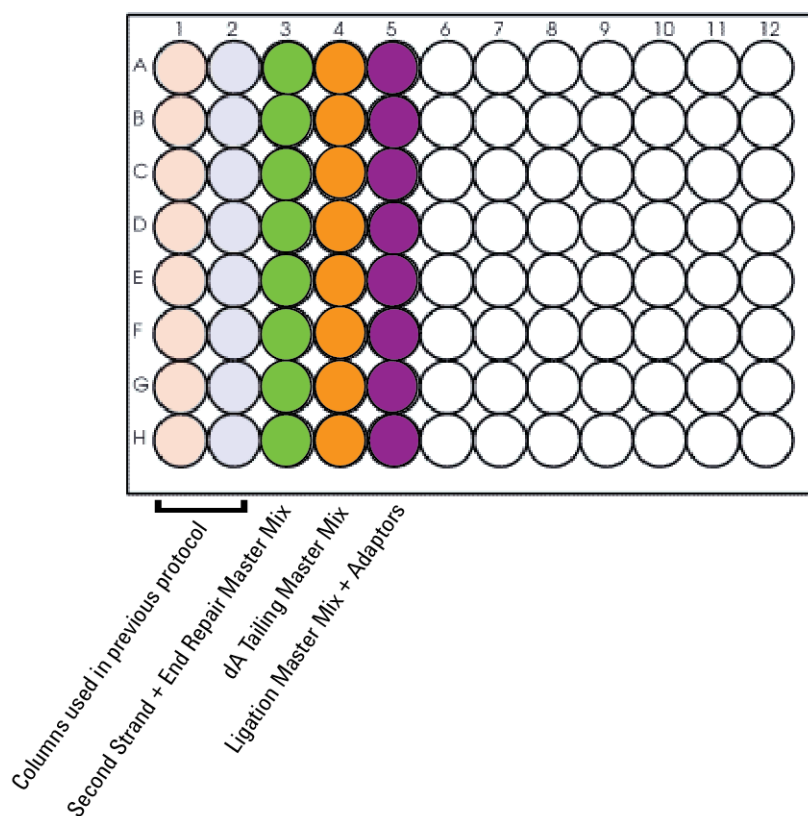


Figure 6 Configuration of the master mix source plate for LibraryPrep_RNASeq_ILM_v1.1.rst

NOTE

If you are using a new DeepWell plate for the Library Prep Master Mix source plate, leave columns 1 and 2 empty and add the PCR Master Mix to columns 3 to 5 of the new plate.

13 Seal the master mix source plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 1.0 sec.

3 Sample Preparation

Step 3. Prepare cDNA libraries for Hybridization

14 Centrifuge the plate for 30 seconds to drive the well contents off the walls and plate seal and to eliminate any bubbles. Keep the master mix source plate on ice.

NOTE

The presence of bubbles in source plate solutions may cause inaccurate volume transfer by the Bravo liquid handling platform. Ensure that the source plate is sealed and centrifuged prior to use in a run.

Load the Agilent NGS Workstation

15 Load the Labware MiniHub according to [Table 30](#), using the plate orientations shown in [Figure 4](#).

Table 30 Initial MiniHub configuration for LibraryPrep_RNASeq_ILM_v1.1.rst

Vertical Shelf Position	Cassette 1	Cassette 2	Cassette 3	Cassette 4
Shelf 5 (Top)	Empty	Empty Nunc DeepWell plate	Empty Nunc DeepWell plate	Empty
Shelf 4	Empty	Empty Eppendorf twin.tec plate	Empty Eppendorf twin.tec plate	Empty
Shelf 3	Empty	Empty	Empty Eppendorf twin.tec plate	Empty
Shelf 2	Waste tip box (retained from mRNA_Purification protocol)	Nuclease-free water reservoir from step 7	AMPure XP beads in Nunc DeepWell plate from step 6	Empty
Shelf 1 (Bottom)	Clean tip box (retained from mRNA_Purification protocol)	70% ethanol reservoir from step 8	Empty	Empty tip box

16 Load the Bravo deck according to [Table 31](#).

Table 31 Initial Bravo deck configuration for LibraryPrep_RNASeq_ILM_v1.1.rst

Location	Content
1	Empty waste reservoir (Axygen 96 Deep Well Plate, square wells)
6	Empty Eppendorf twin.tec plate, oriented with well A1 in the upper-left
7	Purified first-strand cDNA samples in Eppendorf twin.tec plate, oriented with well A1 in the upper-left
9	Library Prep Master Mix Source Plate (Nunc DeepWell), unsealed and seated on silver insert

17 Load the BenchCel Microplate Handling Workstation according to [Table 32](#).

Table 32 Initial BenchCel configuration for LibraryPrep_RNASeq_ILM_v1.1.rst

No. of Columns Processed	Rack 1	Rack 2	Rack 3	Rack 4
1	2 Tip boxes	Empty	Empty	Empty
2	3 Tip boxes	Empty	Empty	Empty
3	4 Tip boxes	Empty	Empty	Empty
4	5 Tip boxes	Empty	Empty	Empty
6	7 Tip boxes	Empty	Empty	Empty
12	11 Tip boxes	3 Tip boxes	Empty	Empty

Run VWorks runset LibraryPrep_RNASeq_ILM_v1.1.rst

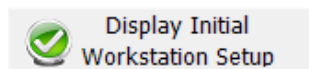
18 On the SureSelect setup form, under **Select Protocol to Run**, select **LibraryPrep_RNASeq_ILM_v1.1.rst**.

19 Select the number of columns of samples to be processed. Runs must include 1, 2, 3, 4, 6, or 12 columns.

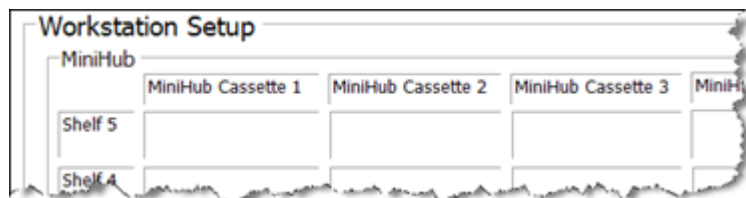
3 Sample Preparation

Step 3. Prepare cDNA libraries for Hybridization

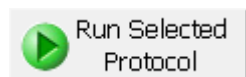
20 Click **Display Initial Workstation Setup**.



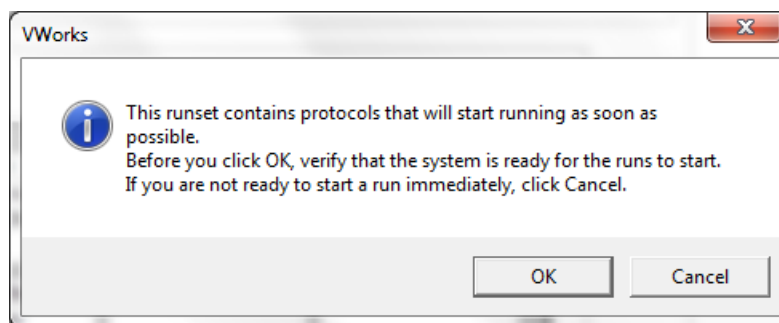
21 Verify that the NGS workstation has been set up as displayed in the Workstation Setup region of the form.



22 When verification is complete, click **Run Selected Protocol**.



23 When ready to begin the run, click **OK** in the following window.



Running the LibraryPrep_RNASeq_ILM_v1.1.rst runset takes approximately 3 hours. Once complete, the purified, adaptor-ligated cDNA samples are located in the Eppendorf plate at position 7 of the Bravo deck.

Stopping Point If you do not continue to the next step, seal the plate and store at -20°C .

Step 4. Amplify cDNA libraries by PCR

In this step, the Agilent NGS Workstation completes the liquid handling steps for PCR amplification of the adaptor-ligated cDNA samples. After the reactions are set up by the workstation, you transfer the PCR plate to a thermal cycler for amplification. The amplification cycle number is based on the initial amount of total RNA sample used for library preparation.

Before you begin, thaw and mix the reagents listed in [Table 33](#) below and keep on ice.

Table 33 Reagents for use in pre-capture PCR protocol

Storage Location	Kit Component	Where Used in Protocol
SureSelect RNA Library Prep, ILM (Pre PCR)*, -20°C	PCR Master Mix [†] (bottle)	page 68
	Uracil DNA Glycosylase (UDG) (tube with yellow cap)	page 68
	SureSelect Primer (tube with brown cap)	page 68
	ILM Reverse PCR Primer [‡] (tube with black cap)	page 68

* May also be labeled as *SureSelect Strand Specific RNA Library Prep, ILM, Box 1*.

† May also be labeled as *RNA Seq PCR Master Mix*.

‡ May also be labeled as *RNA Seq ILM Reverse PCR Primer*.

Prepare the workstation

- 1 Turn on the ThermoCube, set to 0°C, at position 9 of the Bravo deck. Be sure that the chiller reservoir contains at least 300 mL of 25% ethanol.
- 2 Leave tip boxes on shelves 1 and 2 in cassette 1 of the Labware MiniHub from the previous LibraryPrep_RNASeq_v1.0.rst run. Otherwise, clear the remaining positions of the MiniHub and BenchCel of plates and tip boxes.
- 3 Pre-set the temperature of Bravo deck position 6 to 4°C using the Inheco Multi TEC control touchscreen, as described in [Setting the Temperature of Bravo Deck Heat Blocks](#). Bravo deck position 6 corresponds to CPAC 2, position 2 on the Multi TEC control touchscreen.

3 Sample Preparation

Step 4. Amplify cDNA libraries by PCR

Prepare the PCR reaction mix and the master mix source plate

CAUTION

The PCR Master Mix used at this step is highly viscous and thorough mixing prior to use is critical for optimal kit performance. Mix by vortexing at high speed for 5 seconds before removing an aliquot for use and after combining with other solutions as directed in [Table 34](#). Pipetting up and down is not sufficient to mix this reagent.

- 4 Vortex the thawed vial of PCR Master Mix 5 seconds at high speed to ensure homogeneity.
- 5 Prepare the appropriate volume of Pre-capture PCR Reaction Mix, according to [Table 34](#). Mix well by vortexing at high speed and keep on ice.

Table 34 Preparation of Pre-capture PCR Reaction Mix

SureSelect Reagent	Volume for 1 Library	Volume for 1 Column	Volume for 2 Columns	Volume for 3 Columns	Volume for 4 Columns	Volume for 6 Columns	Volume for 12 Columns
Nuclease-free water	5.0 µL	61.5 µL	102.5 µL	143.5 µL	184.5 µL	266.5 µL	533.0 µL
PCR Master Mix	25.0 µL	307.5 µL	512.5 µL	717.5 µL	922.5 µL	1332.5 µL	2665 µL
Uracil DNA Glycosylase (UDG)	1.0 µL	12.3 µL	20.5 µL	28.7 µL	36.9 µL	53.3 µL	106.6 µL
SureSelect Primer (Forward primer)	1.0 µL	12.3 µL	20.5 µL	28.7 µL	36.9 µL	53.3 µL	106.6 µL
ILM Reverse PCR Primer	1.0 µL	12.3 µL	20.5 µL	28.7 µL	36.9 µL	53.3 µL	106.6 µL
Total Volume	33 µL	405.9 µL	676.5 µL	947.1 µL	1217.7 µL	1758.9 µL	3517.8 µL

- 6 Using the same Nunc DeepWell master mix source plate that was used for the LibraryPrep_RNASeq_ILM_v1.1.rst run, add the volume of PCR Reaction Mix indicated in Table 35 to all wells of column 6 of the master mix source plate. The final configuration of the master mix source plate is shown in Figure 7.

Table 35 Preparation of the Master Mix Source Plate for Pre-CapturePCR_RNASeq_ILM_v1.0.pro

Master Mix Solution	Position on Source Plate	Volume of Master Mix added per Well of Nunc Deep Well Source Plate					
		1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
PCR Reaction Mix	Column 6 (A6-H6)	48.6 μ L	80.4 μ L	114.3 μ L	148.1 μ L	215.7 μ L	435.6 μ L

NOTE

If you are using a new DeepWell plate for the pre-capture PCR source plate, leave columns 1 to 5 empty and add the PCR Master Mix to column 6 of the new plate.

3 Sample Preparation

Step 4. Amplify cDNA libraries by PCR

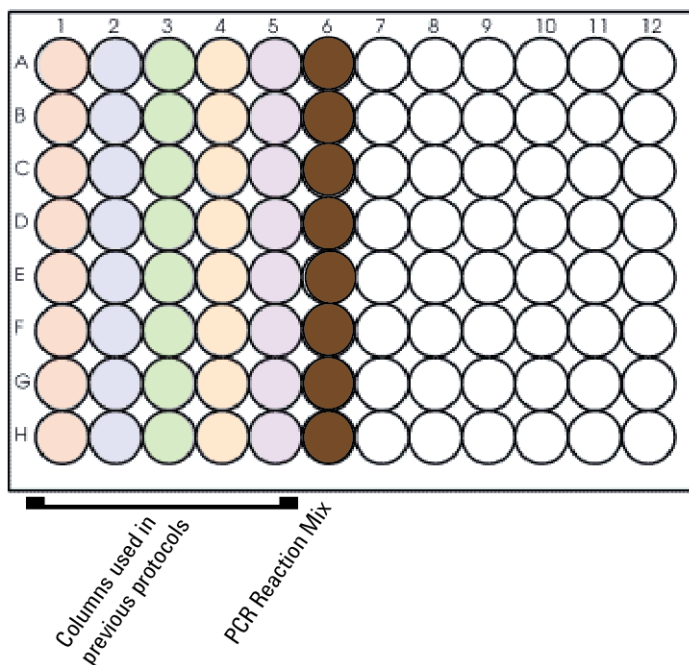


Figure 7 Configuration of the master mix source plate for Pre-CapturePCR_RNASeq_ILM_v1.0.pro. Columns 1-5 were used to dispense master mixes during previous protocols.

- 7 Seal the master mix source plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 1.0 sec.
- 8 Centrifuge the plate for 30 seconds to drive the well contents off the walls and plate seal and to eliminate any bubbles.

NOTE

The presence of bubbles in source plate solutions may cause inaccurate volume transfer by the Bravo liquid handling platform. Ensure that the source plate is sealed and centrifuged prior to use in a run.

Load the Agilent NGS Workstation

9 Load the Labware MiniHub according to [Table 36](#), using the plate orientations shown in [Figure 4](#).

Table 36 Initial MiniHub configuration for Pre-CapturePCR_RNASeq_ILM_v1.0.pro

Vertical Shelf Position	Cassette 1	Cassette 2	Cassette 3	Cassette 4
Shelf 5 (Top)	Empty	Empty	Empty	Empty
Shelf 4	Empty	Empty	Empty	Empty
Shelf 3	Empty	Empty	Empty	Empty
Shelf 2	Waste tip box (retained from Library Prep protocol)	Empty	Empty	Empty
Shelf 1 (Bottom)	Clean tip box (retained from Library Prep protocol)	Empty	Empty	Empty tip box

10 Load the Bravo deck according to [Table 37](#).

Table 37 Initial Bravo deck configuration for Pre-CapturePCR_RNASeq_ILM_v1.0.pro

Location	Content
6	Empty PCR plate seated on red insert (PCR plate type must be specified on setup form under step 2)
7	Prepped cDNA samples in Eppendorf twin.tec plate, oriented with well A1 in the upper-left
9	Master mix source plate (Nunc DeepWell), unsealed and seated on silver insert

3 Sample Preparation

Step 4. Amplify cDNA libraries by PCR

11 Load the BenchCel Microplate Handling Workstation according to [Table 38](#).

Table 38 Initial BenchCel configuration for Pre-CapturePCR_RNASeq_ILM_v1.0.pro

No. of Columns Processed	Rack 1	Rack 2	Rack 3	Rack 4
1	1 Tip box	Empty	Empty	Empty
2	1 Tip box	Empty	Empty	Empty
3	1 Tip box	Empty	Empty	Empty
4	1 Tip box	Empty	Empty	Empty
6	1 Tip box	Empty	Empty	Empty
12	1 Tip box	Empty	Empty	Empty

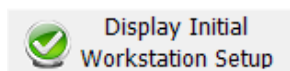
Run VWorks protocol Pre-CapturePCR_RNASeq_ILM_v1.0.pro

12 On the SureSelect setup form, under **Select Protocol to Run**, select **Pre-CapturePCR_RNASeq_ILM_v1.0.pro**.

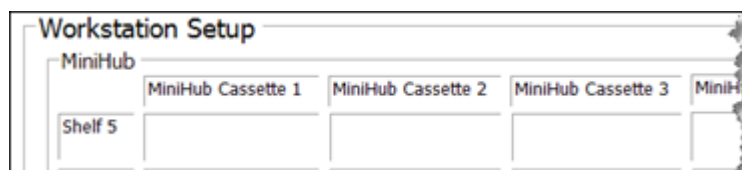
13 Under **Select PCR plate labware for Thermal Cycling**, select the specific type of PCR plate that was loaded on Bravo deck position 6. The plate type selected must be compatible with the thermal cycler to be used for amplification.

14 Select the number of columns of samples to be processed. Runs must include 1, 2, 3, 4, 6, or 12 columns.

15 Click **Display Initial Workstation Setup**.



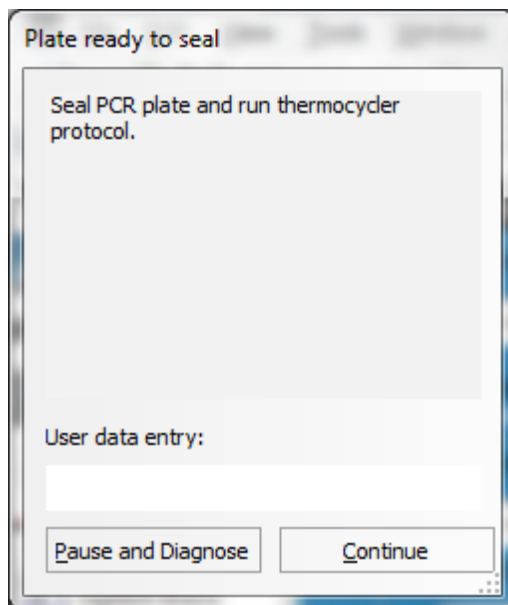
16 Verify that the NGS workstation has been set up as displayed in the Workstation Setup region of the form.



17 When verification is complete, click **Run Selected Protocol**.



18 Running the Pre-CapturePCR_RNASeq_ILM_v1.0.pro protocol takes approximately 15 minutes. Once complete, the PCR-ready samples, containing prepped DNA and PCR master mix are located in the PCR plate at position 6 of the Bravo deck and you will see the following prompt:



- a Remove the PCR plate from position 6 of the Bravo deck and seal the plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 3.0 seconds.
- b Centrifuge the plate for 30 seconds to drive the well contents off the walls and plate seal and to eliminate air bubbles.
- c Transfer the plate to the thermal cycler (with the heated lid ON) and run the program in [Table 39](#).

3 Sample Preparation

Step 4. Amplify cDNA libraries by PCR

Table 39 Thermal cycler program for mRNA Library PCR indexing

Segment	Number of Cycles	Temperature	Time
1	1	37°C	15 minutes
2	1	95°C	2 minutes
3	9–13 cycles (see Table 40)	95°C	30 seconds
		65°C	30 seconds
		72°C	1 minute
4	1	72°C	5 minutes
5	1	4°C	Hold

Table 40 mRNA Library PCR indexing cycle number recommendations

Amount of total RNA used for library prep	Cycle Number
200 ng–2 µg	11–13
2.1 µg–4 µg	9–11

Step 5. Purify amplified DNA using AMPure XP beads

In this step, the Agilent NGS Workstation transfers AMPure XP beads and amplified cDNA libraries to a Nunc DeepWell plate and then collects and washes the bead-bound DNA.

Prepare the workstation and reagents

- 1 Clear the Labware MiniHub and BenchCel of all plates and tip boxes.
- 2 Gently wipe down the Labware MiniHub, Bravo deck, and BenchCel with a DNA Away decontamination wipe.
- 3 Turn on the ThermoCube, set to 0°C, at position 9 of the Bravo deck. Be sure that the chiller reservoir contains at least 300 mL of 25% ethanol.
- 4 Let the AMPure XP beads come to room temperature for at least 30 minutes. *Do not freeze the beads at any time.*
- 5 Mix the bead suspension well so that the reagent appears homogeneous and consistent in color.
- 6 Prepare a Nunc DeepWell source plate for the beads by adding 65 µL of homogeneous AMPure XP beads per well, for each well to be processed.
- 7 Prepare a Thermo Scientific reservoir containing 15 mL of nuclease-free water.
- 8 Prepare a separate Thermo Scientific reservoir containing 45 mL of freshly-prepared 70% ethanol.

3 Sample Preparation

Step 5. Purify amplified DNA using AMPure XP beads

9 Load the Labware MiniHub according to [Table 41](#), using the plate orientations shown in [Figure 4](#).

Table 41 Initial MiniHub configuration for AMPureXP_v1.1.pro:Pre-Capture PCR

Vertical Shelf Position	Cassette 1	Cassette 2	Cassette 3	Cassette 4
Shelf 5 (Top)	Empty Nunc DeepWell plate	Empty	Empty	Empty
Shelf 4	Empty	Empty	Empty	Empty
Shelf 3	Empty	Empty Eppendorf twin.tec Plate	Empty	Empty
Shelf 2	Empty	Nuclease-free water reservoir from step 7	AMPure XP beads in Nunc DeepWell plate from step 6	Empty
Shelf 1 (Bottom)	Empty	70% ethanol reservoir from step 8	Empty	Empty tip box

10 Load the Bravo deck according to [Table 42](#).

Table 42 Initial Bravo deck configuration for AMPureXP_v1.1.pro:Pre-Capture PCR

Location	Content
1	Empty waste reservoir (Axygen 96 Deep Well Plate, square wells)
9	Amplified cDNA samples in PCR plate, unsealed and seated on red insert (PCR plate type must be specified on setup form under step 2)

Step 5. Purify amplified DNA using AMPure XP beads

11 Load the BenchCel Microplate Handling Workstation according to Table 43.

Table 43 Initial BenchCel configuration for AMPureXP_v1.1.pro:Pre-Capture PCR

No. of Columns Processed	Rack 1	Rack 2	Rack 3	Rack 4
1	1 Tip box	Empty	Empty	Empty
2	1 Tip box	Empty	Empty	Empty
3	2 Tip boxes	Empty	Empty	Empty
4	2 Tip boxes	Empty	Empty	Empty
6	3 Tip boxes	Empty	Empty	Empty
12	6 Tip boxes	Empty	Empty	Empty

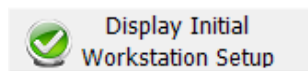
Run VWorks protocol AMPureXP_v1.1.pro:Pre-Capture PCR

12 On the SureSelect setup form, under **Select Protocol to Run**, select **AMPureXP_v1.1.pro:Pre-Capture PCR**.

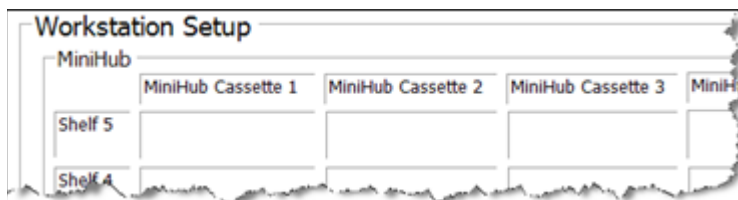
13 Under **Select PCR plate labware for Thermal Cycling**, select the specific type of PCR plate that was loaded on Bravo deck position 9.

14 Select the number of columns of samples to be processed. Runs must include 1, 2, 3, 4, 6, or 12 columns.

15 Click **Display Initial Workstation Setup**.



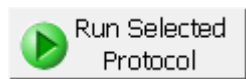
16 Verify that the NGS workstation has been set up as displayed in the Workstation Setup region of the form.



3 Sample Preparation

Step 5. Purify amplified DNA using AMPure XP beads

17 When verification is complete, click **Run Selected Protocol**.



The purification protocol takes approximately 45 minutes. When complete, the purified DNA samples are in the Eppendorf plate located on Bravo deck position 7.

Stopping Point If you do not continue to the next step, seal the plate and store at -20°C .

Step 6. Assess library DNA quantity and quality

Option 1: Analysis using the Agilent 2100 Bioanalyzer and DNA 1000 Assay

Use a Bioanalyzer DNA 1000 chip and reagent kit. Perform the assay according to the [Agilent DNA 1000 Kit Guide](#).

- 1 Set up the 2100 Bioanalyzer as instructed in the reagent kit guide.
- 2 Seal the sample plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 1.0 sec.
- 3 Vortex the plate to mix samples in each well, then centrifuge the plate for 30 seconds to drive the well contents off the walls and plate seal.
- 4 Prepare the chip, samples and ladder as instructed in the reagent kit guide, using 1 μL of each sample for the analysis.
- 5 Load the prepared chip into the 2100 Bioanalyzer and start the run within five minutes after preparation.
- 6 Measure the concentration of the library (ng/ μL) by integrating under the peak at approximately 180 to 550 bp. For accurate quantification, make sure that the concentration falls within the linear range of the assay.

A sample electropherogram is shown in [Figure 8](#).

Stopping Point

If you do not continue to the next step, seal the plate and store at -20°C .

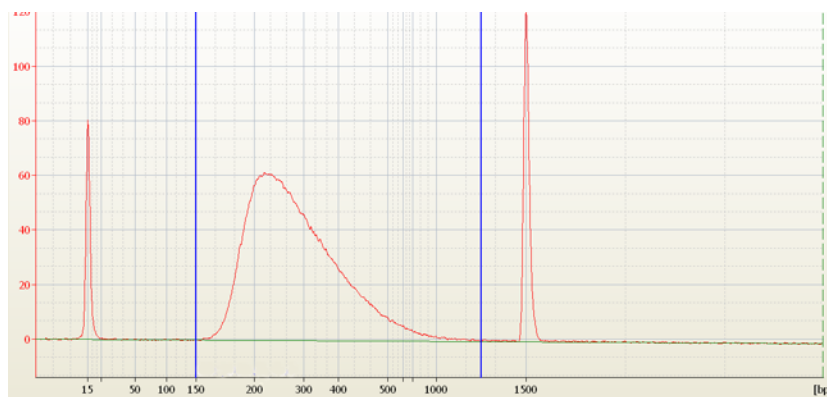


Figure 8 Analysis of amplified library DNA using a DNA 1000 assay.

3 Sample Preparation

Step 6. Assess library DNA quantity and quality

Option 2: Analysis using Agilent 4200 TapeStation and D1000 ScreenTape

Use a D1000 ScreenTape and associated reagent kit to analyze the amplified libraries. For more information to do this step, see the [Agilent D1000 Assay Quick Guide](#).

- 1 Seal the DNA sample plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 1.0 sec.
- 2 Vortex the plate to mix samples in each well, then centrifuge the plate for 30 seconds to drive the well contents off the walls and plate seal.
- 3 Prepare the TapeStation samples as instructed in the instrument user manual. Use 1 µL of each amplified library DNA sample diluted with 3 µL of D1000 sample buffer for the analysis.

CAUTION

For accurate quantitation, make sure to thoroughly mix the combined DNA and sample buffer by vortexing the assay plate or tube strip for 1 minute on the IKA MS3 vortex mixer provided with the 4200 TapeStation system before loading the samples.

If no IKA MS3 vortex mixer is available, ensure that the samples are mixed thoroughly by vortexing for 10 seconds at high speed on a manual vortex mixer.

-
- 4 Load the sample plate or tube strips from [step 3](#), the D1000 ScreenTape, and loading tips into the TapeStation as instructed in the [Assay Quick Guide](#). Start the run.
 - 5 For each sample, measure the concentration of the library (ng/µL) by integrating under the peak at approximately 180 to 550 bp. A sample electropherogram is shown in [Figure 9](#).

Stopping Point If you do not continue to the next step, seal the plate and store at -20°C.

Step 6. Assess library DNA quantity and quality

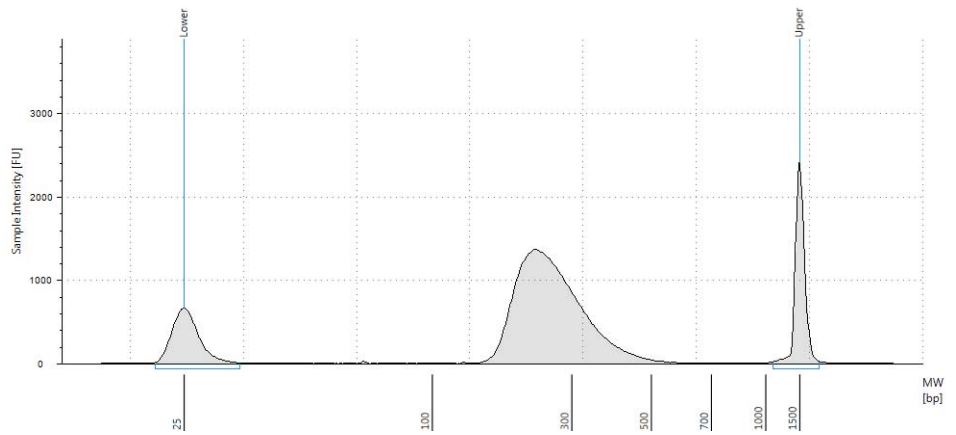
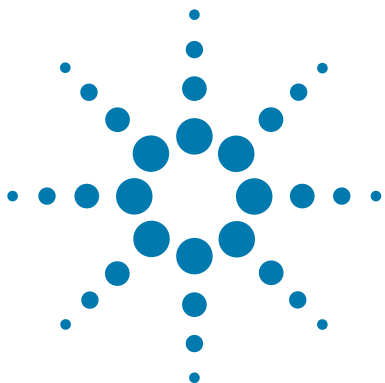


Figure 9 Analysis of amplified library DNA using the D1000 ScreenTape.

3 Sample Preparation

Step 6. Assess library DNA quantity and quality



4 Hybridization

Step 1. Aliquot prepped DNA libraries for hybridization 84

Step 2. Hybridize the DNA library to the Probe 87

Step 3. Capture the hybridized DNA 100

This chapter describes the steps to combine the prepped library with the blocking agents and the SureSelect or ClearSeq RNA Target Enrichment Probe. Each cDNA library sample must be hybridized and captured individually prior to addition of the indexing tag by PCR.

CAUTION

The ratio of probe to prepped library is critical for successful capture.

CAUTION

You must avoid evaporation from the small volumes of the capture during the 24 hour incubation.

If you want to use a duration of hybridization >24 hours, first test the conditions. Incubate 35 μ L of SureSelect Hybridization Buffer (without DNA) at 65°C for 24 hours (or longer, if applicable) as a test. Include buffer in each well that you might use, including those in the center and those on the edges. Check that you do not get extensive evaporation. Evaporation should not exceed 3 to 4 μ L.



4 Hybridization

Step 1. Aliquot prepped DNA libraries for hybridization

Step 1. Aliquot prepped DNA libraries for hybridization

For each sample library prepared, do one hybridization and capture. Do not pool samples at this stage.

Each hybridization reaction will contain 100 ng of the appropriate prepped cDNA sample. Before starting the hybridization step, you must create a table containing instructions for the Agilent NGS Workstation indicating the volume of each sample required for a 100-ng aliquot.

- 1 Create a .csv (comma separated value) file with the headers shown in [Figure 10](#). The header text must not contain spaces. The table may be created using a spreadsheet application, such as Microsoft Excel software, and then saved in .csv format. The file must include rows for all 96 wells of the plate.
- 2 Enter the information requested in the header for each DNA sample.
 - In the SourceBC field, enter the sample plate description or barcode. The SourceBC field contents must be identical for all rows.
 - In the SourceWell and DestinationWell fields, enter each well position for the plate. SourceWell and DestinationWell field contents must be identical for a given sample.
 - In the Volume field, enter the volume (in μL) equivalent to 100 ng DNA for each sample. These values are determined from the concentration values obtained from Bioanalyzer or TapeStation traces in the previous section. For all empty wells on the plate, enter the value 0, as shown in [Figure 10](#); **do not delete rows for empty wells**.

	A	B	C	D
1	SourceBC	SourceWell	DestinationWell	Volume
2	SamplePlateXYZ	A1	A1	5.35
3	SamplePlateXYZ	B1	B1	4.28
4	SamplePlateXYZ	C1	C1	4.76
5	SamplePlateXYZ	D1	D1	5.19
6	SamplePlateXYZ	E1	E1	5.49
7	SamplePlateXYZ	F1	F1	4.86
8	SamplePlateXYZ	G1	G1	5.05
9	SamplePlateXYZ	H1	H1	4.37
10	SamplePlateXYZ	A2	A2	0
11	SamplePlateXYZ	B2	B2	0
12	SamplePlateXYZ	C2	C2	0
13	SamplePlateXYZ	D2	D2	0

Figure 10 Sample spreadsheet for 100-ng sample aliquot for 1-column run.

NOTE

You can find a sample spreadsheet in the directory **C: > VWorks Workspace > NGS Option B > XT_RNA_ILM > Aliquot Library Input Files > 100ng_transfer_full_plate_template.xlsx**.

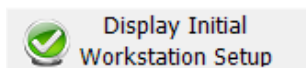
The 100ng_transfer_full_plate_template.xlsx file may be copied and used as a template for creating the .csv files for each Aliquot_Libraries_v1.0.pro run. If you are using the sample file as a template for runs with fewer than 12 columns, be sure to retain rows for all 96 wells, and populate the Volume column with 0 for unused wells.

- 3 Load the .csv file onto the PC containing the VWorks software into a suitable folder, such as **C: > VWorks Workspace > NGS Option B > XT_RNA_ILM > Aliquot Library Input Files**.
- 4 Turn on the chiller, set to 0°C, at position 9 of the Bravo deck. Be sure that the chiller reservoir contains at least 300 mL of 25% ethanol.
- 5 Load the Bravo deck according to [Table 44](#).

Table 44 Initial Bravo deck configuration for Aliquot_Libraries_v1.0.pro

Location	Content
5	Empty PCR plate seated on red insert (PCR plate type must be specified on setup form under step 2)
6	Empty tip box
8	New tip box
9	Prepped library DNA in Eppendorf plate (oriented with well A1 in the upper-left)

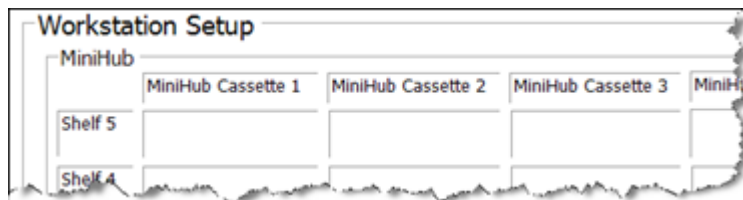
- 6 On the SureSelect setup form, under **Select Protocol to Run**, select **Aliquot_Libraries_v1.0.pro**.
- 7 Click **Display Initial Workstation Setup**.



4 Hybridization

Step 1. Aliquot prepped DNA libraries for hybridization

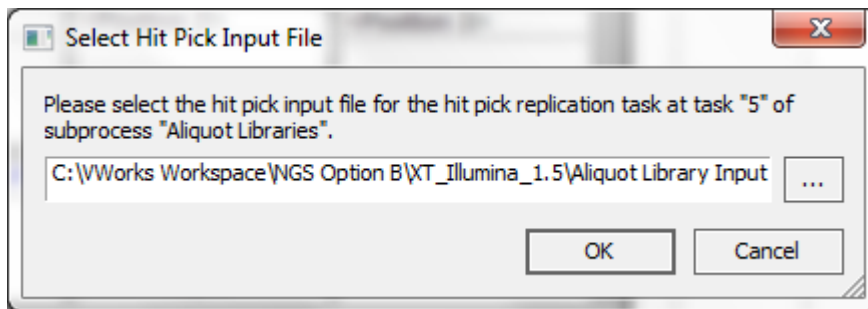
- 8 Verify that the NGS workstation has been set up as displayed in the Workstation Setup region of the form.



- 9 When verification is complete, click **Run Selected Protocol**.



- 10 When prompted by the dialog below, browse to the .csv file created for the source plate of the current run, and then click **OK** to start the run.



The library aliquoting protocol takes approximately 1 hour for 96 samples. When complete, the 100-ng samples are in the PCR plate located on Bravo deck position 5.

- 11 Remove the 100-ng sample plate from the Bravo deck and use a vacuum concentrator to dry the sample at $\leq 45^{\circ}\text{C}$.
- 12 Reconstitute each dried sample with 3.4 μL of nuclease-free water to bring the final concentration to 29.4 ng/ μL . Pipette up and down along the sides of each well for optimal recovery.
- 13 Seal the plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 1.0 sec.
- 14 Vortex the plate for 30 seconds to ensure complete reconstitution, then centrifuge the plate for 1 minute to drive the well contents off the walls and plate seal.

Step 2. Hybridize the DNA library to the Probe

In this step, the Agilent NGS Workstation completes the liquid handling steps to prepare for hybridization. Afterward, you transfer the sample plate to a thermal cycler, held at 65°C, to allow hybridization of the prepared cDNA samples to one or more RNA Target Enrichment Probes.

Prepare the workstation

- 1 Clear the Labware MiniHub and BenchCel of all plates and tip boxes.
- 2 Gently wipe down the Labware MiniHub, Bravo decks, and BenchCel with a DNA Away decontamination wipe.
- 3 Turn on the ThermoCube and set to 25°C for position 9 of the Bravo deck.
- 4 Place the silver Nunc DeepWell plate insert on position 9 of the Bravo deck. This insert is required to facilitate heat transfer to DeepWell source plate wells during the Hybridization protocol.

4 Hybridization

Step 2. Hybridize the DNA library to the Probe

Prepare the SureSelect Block master mix

- 5 Prepare the appropriate volume of SureSelect Block master mix, on ice, as indicated in [Table 45](#).

Table 45 Preparation of SureSelect Block Master Mix

SureSelect ^{XT} Reagent	Volume for 1 Library	Volume for 1 Column	Volume for 2 Columns	Volume for 3 Columns	Volume for 4 Columns	Volume for 6 Columns	Volume for 12 Columns
Nuclease-free water	6.0 µL	76.5 µL	127.5 µL	178.5 µL	229.5 µL	331.5 µL	663.0 µL
SureSelect Indexing Block 1 (green cap)	2.5 µL	31.9 µL	53.1 µL	74.4 µL	95.6 µL	138.1 µL	276.3 µL
SureSelect Block 2 (blue cap)	2.5 µL	31.9 µL	53.1 µL	74.4 µL	95.6 µL	138.1 µL	276.3 µL
SureSelect Indexing Block 3 (brown cap)	0.6 µL	7.7 µL	12.8 µL	17.9 µL	23.0 µL	33.2 µL	66.3 µL
Total Volume	11.6 µL	147.9 µL	246.5 µL	345.1 µL	443.7 µL	640.9 µL	1281.9 µL

Prepare one or more Capture Library Master Mixes

- 6 Prepare the appropriate volume of Capture Library Master Mix for each of the probes that will be used for hybridization as indicated in [Table 46](#) to [Table 49](#). Mix the components by pipetting. Keep the master mixes on ice during preparation and aliquoting.

NOTE

Each row of the prepped cDNA sample plate may be hybridized to a different Probe.

For runs that use a single Probe for all rows of the plate, prepare a single Capture Library Master Mix as described in Step a ([Table 46](#) or [Table 47](#)) below.

For runs that use different Probes for individual rows, prepare each of the multiple Capture Library Master Mixes as described in Step b ([Table 48](#) or [Table 49](#)) below.

Step 2. Hybridize the DNA library to the Probe

- a For runs that use a single Probe for all rows**, prepare the Capture Library Master Mix as listed in [Table 46](#) or [Table 47](#), based on the Mb target size of your design.

Table 46 Preparation of Capture Library Master Mix for target sizes <3.0 Mb, 8 rows of wells

Target size <3.0 Mb							
SureSelect ^{XT} Reagent	Volume for 1 Library	Volume for 1 Column	Volume for 2 Columns	Volume for 3 Columns	Volume for 4 Columns	Volume for 6 Columns	Volume for 12 Columns
Nuclease-free water	4.5 µL	76.5 µL	114.8 µL	153.0 µL	191.3 µL	306.0 µL	612.0 µL
RNase Block (purple cap)	0.5 µL	8.5 µL	12.8 µL	17.0 µL	21.3 µL	34.0 µL	68.0 µL
SureSelect or ClearSeq Probe	2.0 µL	34.0 µL	51.0 µL	68.0 µL	85.0 µL	136.0 µL	272.0 µL
Total Volume	7.0 µL	119.0 µL	178.6 µL	238.0 µL	297.6 µL	476.0 µL	852.0 µL

Table 47 Preparation of Capture Library Master Mix for target sizes >3.0 Mb, 8 rows of wells

Target size >3.0 Mb							
SureSelect ^{XT} Reagent	Volume for 1 Library	Volume for 1 Column	Volume for 2 Columns	Volume for 3 Columns	Volume for 4 Columns	Volume for 6 Columns	Volume for 12 Columns
Nuclease-free water	1.5 µL	25.5 µL	38.3 µL	51.0 µL	63.8 µL	102.0 µL	204.0 µL
RNase Block (purple cap)	0.5 µL	8.5 µL	12.8 µL	17.0 µL	21.3 µL	34.0 µL	68.0 µL
SureSelect or ClearSeq Probe	5.0 µL	85.0 µL	127.5 µL	170.0 µL	212.5 µL	340.0 µL	680.0 µL
Total Volume	7.0 µL	119.0 µL	178.6 µL	238.0 µL	297.6 µL	476.0 µL	852.0 µL

4 Hybridization

Step 2. Hybridize the DNA library to the Probe

- b For runs that use different Probes in individual rows**, prepare each Capture Library Master Mix as listed in [Table 48](#) or [Table 49](#), based on the Mb target size of each design. The volumes listed in [Table 48](#) and [Table 49](#) are for a single row of sample wells. If a given Probe will be hybridized in multiple rows, multiply each of the values below by the number of rows assigned to that Probe.

Table 48 Preparation of Capture Library Master Mix for target sizes <3.0 Mb, single row of wells

Target size <3.0 Mb							
SureSelect ^{XT} Reagent	Volume for 1 Library	Volume for 1 Column	Volume for 2 Columns	Volume for 3 Columns	Volume for 4 Columns	Volume for 6 Columns	Volume for 12 Columns
Nuclease-free water	4.5 µL	9.0 µL	13.8 µL	18.6 µL	23.3 µL	37.7 µL	75.9 µL
RNase Block (purple cap)	0.5 µL	1.0 µL	1.5 µL	2.1 µL	2.6 µL	4.2 µL	8.4 µL
SureSelect or ClearSeq Probe	2.0 µL	4.0 µL	6.1 µL	8.3 µL	10.4 µL	16.8 µL	33.8 µL
Total Volume	7.0 µL	14.0 µL	21.4 µL	28.9 µL	36.3 µL	58.6 µL	118.1 µL

Table 49 Preparation of Capture Library Master Mix for target sizes >3.0 Mb, single row of wells

Target size >3.0 Mb							
SureSelect ^{XT} Reagent	Volume for 1 Library	Volume for 1 Column	Volume for 2 Columns	Volume for 3 Columns	Volume for 4 Columns	Volume for 6 Columns	Volume for 12 Columns
Nuclease-free water	1.5 µL	3.0 µL	4.6 µL	6.2 µL	7.8 µL	12.6 µL	25.3 µL
RNase Block (purple cap)	0.5 µL	1.0 µL	1.5 µL	2.1 µL	2.6 µL	4.2 µL	8.4 µL
SureSelect or ClearSeq Probe	5.0 µL	10.0 µL	15.3 µL	20.6 µL	25.9 µL	41.9 µL	84.4 µL
Total Volume	7.0 µL	14.0 µL	21.4 µL	28.9 µL	36.3 µL	58.6 µL	118.1 µL

Prepare the Hybridization Buffer master mix

- 7 Prepare the appropriate volume of Hybridization Buffer Master Mix, at room temperature, as indicated in [Table 50](#).

Table 50 Preparation of Hybridization Buffer Master Mix

SureSelect ^{XT} Reagent	Volume for 1 Column	Volume for 2 Columns	Volume for 3 Columns	Volume for 4 Columns	Volume for 6 Columns	Volume for 12 Columns
SureSelect Hyb 1 (bottle)	140.9 µL	197.3 µL	250.0 µL	310.1 µL	422.8 µL	789.3 µL
SureSelect Hyb 2 (red cap)	5.6 µL	7.9 µL	10.0 µL	12.4 µL	16.9 µL	31.6 µL
SureSelect Hyb 3 (yellow cap)	56.4 µL	78.9 µL	100.0 µL	124.0 µL	169.1 µL	315.7 µL
SureSelect Hyb 4 (black cap)	73.3 µL	102.6 µL	130.0 µL	161.2 µL	219.9 µL	410.4 µL
Total Volume	276.2 µL	386.7 µL	490.0 µL	607.7 µL	828.7 µL	1547 µL

- 8 If precipitate forms, warm the hybridization buffer at 65°C for 5 minutes.

4 Hybridization

Step 2. Hybridize the DNA library to the Probe

Prepare the master mix source plate

- 9 In a Nunc DeepWell plate, prepare the master mix source plate containing the master mixes prepared in [step 5](#) to [step 7](#) at room temperature. Add the volumes indicated in [Table 51](#) of each master mix to each well of the indicated column of the Nunc DeepWell plate. When using multiple Capture Libraries in a run, add each Capture Library Master Mix to the appropriate row(s) of the Nunc DeepWell plate. The final configuration of the master mix source plate is shown in [Figure 11](#).

Table 51 Preparation of the Master Mix Source Plate for SureSelectHybridization_v1.0.pro

Master Mix Solution	Position on Source Plate	Volume of Master Mix added per Well of Nunc Deep Well Source Plate					
		1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
Block Master Mix	Column 1 (A1-H1)	17.0 µL	29.4 µL	41.7 µL	54.0 µL	78.7 µL	158.8 µL
Capture Library Master Mix	Column 2 (A2-H2)	14.0 µL	21.4 µL	28.9 µL	36.3 µL	51.2 µL	99.5 µL
Hybridization Buffer Master Mix	Column 3 (A3-H3)	30.5 µL	44.3 µL	57.2 µL	71.9 µL	99.5 µL	189.3 µL

Step 2. Hybridize the DNA library to the Probe

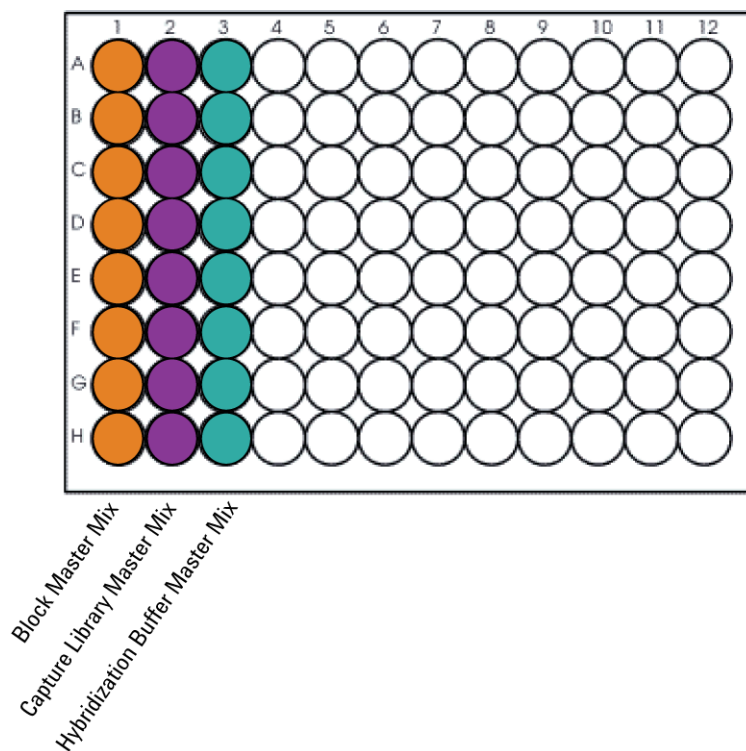


Figure 11 Configuration of the master mix source plate for SureSelectHybridization_v1.0.pro.

- 10** Seal the master mix source plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 1.0 sec.
- 11** Centrifuge the plate for 30 seconds to drive the well contents off the walls and plate seal and to eliminate any bubbles. Keep the master mix plate at room temperature.

4 Hybridization

Step 2. Hybridize the DNA library to the Probe

Load the Agilent NGS Workstation

12 Load the Bravo deck according to [Table 52](#).

Table 52 Initial Bravo deck configuration for SureSelectHybridization_v1.0.pro

Location	Content
5	Empty Eppendorf twin.tec plate
6	100-ng aliquots of prepped DNA libraries in PCR plate seated on red insert (PCR plate type must be specified on setup form under step 2)
8	Empty tip box
9	Hybridization Master Mix source plate seated on silver insert

13 Load the BenchCel Microplate Handling Workstation according to [Table 53](#).

Table 53 Initial BenchCel configuration for SureSelectHybridization_v1.0.pro

No. of Columns Processed	Rack 1	Rack 2	Rack 3	Rack 4
1	1 Tip box	Empty	Empty	Empty
2	1 Tip box	Empty	Empty	Empty
3	2 Tip boxes	Empty	Empty	Empty
4	2 Tip boxes	Empty	Empty	Empty
6	3 Tip boxes	Empty	Empty	Empty
12	5 Tip boxes	Empty	Empty	Empty

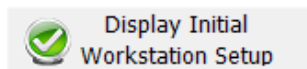
Run VWorks protocol SureSelectHybridization_v1.0.pro

14 On the SureSelect setup form, under **Select Protocol to Run**, select **SureSelectHybridization_v1.0.pro**.

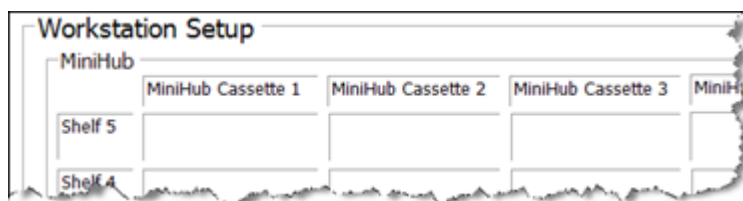
15 Under **Select PCR plate labware for Thermal Cycling**, select the specific type of PCR plate containing the DNA samples at position 6.

16 Select the number of columns of samples to be processed. Runs must include 1, 2, 3, 4, 6, or 12 columns.

17 Click **Display Initial Workstation Setup**.



18 Verify that the NGS workstation has been set up as displayed in the Workstation Setup region of the form.



19 When verification is complete, click **Run Selected Protocol**.

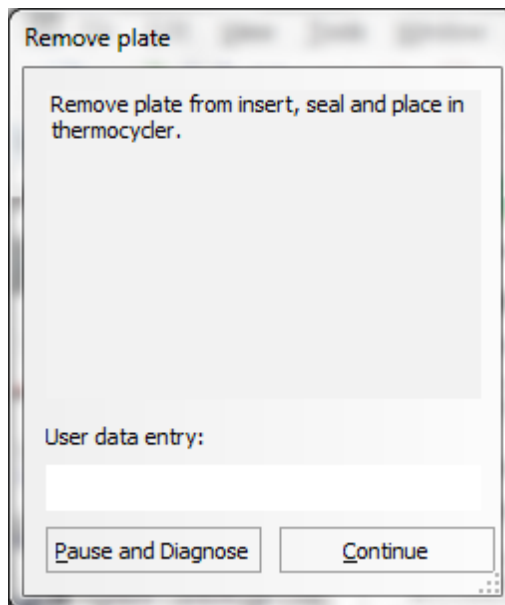


The Agilent NGS Workstation transfers SureSelect Block Master Mix to the prepped DNA-containing wells of the PCR plate. When this process is complete, you will be prompted to transfer the plate to the thermal cycler for sample denaturation prior to hybridization.

4 Hybridization

Step 2. Hybridize the DNA library to the Probe

- 20 When prompted by VWorks as shown below, remove the PCR plate from position 6 of the Bravo deck, leaving the red insert in place.



- 21 Seal the sample plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 3.0 sec.
- 22 Transfer the sealed plate to a thermal cycler and run the following program shown in Table 54. After transferring the plate, click **Continue** on the VWorks screen.

Table 54 Thermal cycler program used for sample denaturation prior to hybridization

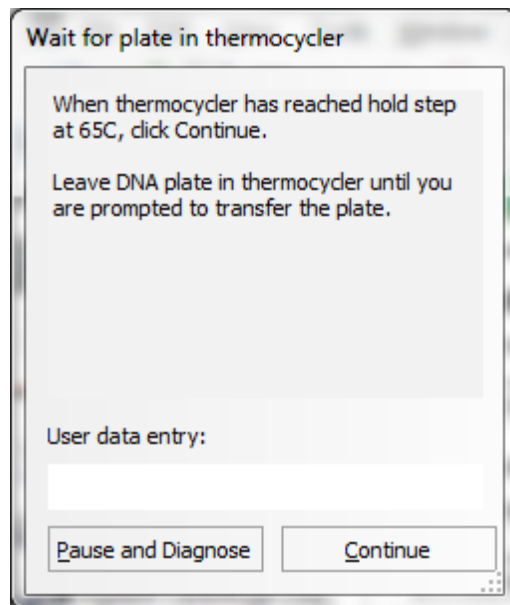
Step	Temperature	Time
Step 1	95°C	5 minutes
Step 2	65°C	Hold

While the sample plate incubates on the thermal cycler, the Agilent NGS Workstation combines aliquots of the Capture Library master mix and Hybridization Buffer master mix.

CAUTION

You must complete [step 23](#) to [step 27](#) quickly, and immediately after being prompted by the VWorks software. It is important that sample temperature remains approximately 65°C during transfers between the Agilent NGS Workstation and thermal cycler.

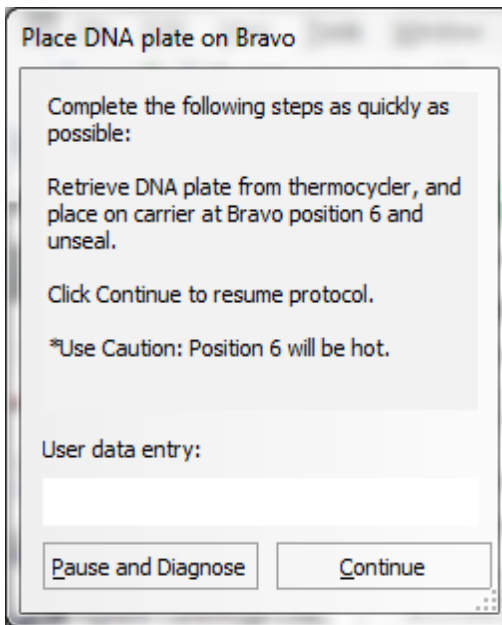
- 23** When the workstation has finished aliquoting the Capture Library and Hybridization Buffer master mixes, you will be prompted by VWorks as shown below. When the thermal cycler reaches the 65°C hold step, click **Continue**. Leave the sample plate in the thermal cycler until you are notified to move it.



4 Hybridization

Step 2. Hybridize the DNA library to the Probe

- 24 When prompted by VWorks as shown below, quickly remove the sample plate from the thermal cycler, unseal the plate carefully to avoid splashing, and transfer the plate to position 6 of the Bravo deck, seated in the red insert. Click **Continue**.



WARNING

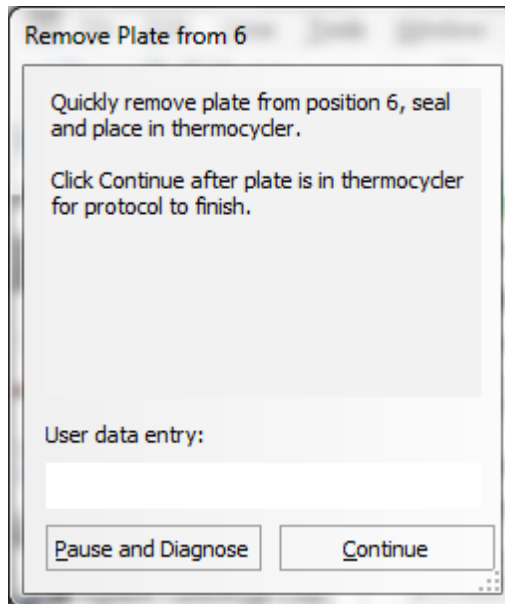
Bravo deck position 6 will be hot.

Use caution when handling components that contact heated deck positions.

The Agilent NGS Workstation transfers the Capture Library-hybridization buffer mixture to the wells of the PCR plate, containing the mixture of prepped DNA samples and blocking agents.

Step 2. Hybridize the DNA library to the Probe

- 25 When prompted by VWorks as shown below, quickly remove the PCR plate from Bravo deck position 6, leaving the insert in place.



- 26 Seal the sample plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 3.0 sec.
- 27 Quickly transfer the plate back to the thermal cycler, held at 65°C. After transferring the plate, click **Continue** on the VWorks screen.
- 28 To finish the VWorks protocol, click **Continue** in the **Unused Tips** and **Empty Tip box** dialogs, and click **Yes** in the **Protocol Complete** dialog.

CAUTION

The temperature of the plate in the thermal cycler should be held at 65°C using a heated lid at 105°C. The lid of the thermal cycler is hot and can cause burns. Use caution when working near the lid.

- 29 Incubate the hybridization mixture in the thermal cycler for 24 hours at 65°C with a heated lid at 105°C. If you are using the SureCycler thermal cycler, place a compression mat over the PCR plate before closing the thermal cycler lid for the 24-hour incubation period.

Step 3. Capture the hybridized DNA

In this step, the cDNA-probe hybrids are captured using streptavidin-coated magnetic beads. This step is run immediately after the 24-hour hybridization period.

This step is automated by the NGS workstation using the SureSelectCapture&Wash_v1.0.rst runset, with a total duration of approximately 3 hours. A workstation operator must be present to complete two actions during the runset, at the time points in the table below. The times provided are approximate; each action is completed in response to a VWorks prompt at the appropriate time in the runset.

Table 55

Operator action	Approximate time after run start
Transfer hybridization reactions from thermal cycler to NGS workstation	<5 minutes
Remove PCR plate from red aluminum insert	5-10 minutes

Prepare the workstation

- 1 Clear the Labware MiniHub and BenchCel of all plates and tip boxes.
- 2 Gently wipe down the Labware MiniHub, Bravo decks, and BenchCel with a DNA Away decontamination wipe.
- 3 Pre-set the temperature of Bravo deck position 4 to 66°C using the Inheco Multi TEC control touchscreen, as described in [Setting the Temperature of Bravo Deck Heat Blocks](#). Bravo deck position 4 corresponds to CPAC 2, position 1 on the Multi TEC control touchscreen.

Prepare the Dynabeads M-270 streptavidin beads**CAUTION**

Use only the recommended Dynabeads M-270 Streptavidin Beads for this automated protocol. Use of other streptavidin bead preparations may adversely affect performance and is not supported by Agilent.

- 4 Vigorously resuspend the Dynabeads M-270 Streptavidin magnetic beads on a vortex mixer. The beads settle during storage.
- 5 Wash the magnetic beads.
 - a In a conical vial, combine the components listed in [Table 56](#). The volumes below include the required overage.

Table 56 Components required for magnetic bead washing procedure

Reagent	Volume for 1 Library	Volume for 1 Column	Volume for 2 Columns	Volume for 3 Columns	Volume for 4 Columns	Volume for 6 Columns	Volume for 12 Columns
M-270 Streptavidin bead suspension	50 µL	425 µL	825 µL	1225 µL	1.65 mL	2.5 mL	5.0 mL
SureSelect Binding Buffer	0.2 mL	1.7 mL	3.3 mL	4.9 mL	6.6 mL	10 mL	20 mL
Total Volume	0.25 mL	2.125 mL	4.125 mL	6.125 mL	8.25 mL	12.5 mL	25 mL

- b Mix the beads on a vortex mixer for 5 seconds.
- c Put the vial into a magnetic device, such as the Dynal magnetic separator.
- d Remove and discard the supernatant.
- e Repeat [step a](#) through [step d](#) for a total of 3 washes. (Retain the beads after each wash and combine with a fresh aliquot of the indicated volume of SureSelect Binding Buffer.)

4 Hybridization

Step 3. Capture the hybridized DNA

- 6 Resuspend the beads in SureSelect Binding buffer, according to [Table 57](#) below.

Table 57 Preparation of magnetic beads for SureSelectCapture&Wash_v1.0.rst

Reagent	Volume for 1 Library	Volume for 1 Column	Volume for 2 Columns	Volume for 3 Columns	Volume for 4 Columns	Volume for 6 Columns	Volume for 12 Columns
SureSelect Binding Buffer	0.2 mL	1.7 mL	3.3 mL	4.9 mL	6.6 mL	10 mL	20 mL

- 7 Prepare a Nunc DeepWell source plate for the washed streptavidin bead suspension. For each well to be processed, add 200 μ L of the homogeneous bead suspension to the Nunc DeepWell plate.
- 8 Place the streptavidin bead source plate at position 5 of the Bravo deck.

Prepare capture and wash solution source plates

- 9 Prepare a Thermo Scientific reservoir containing 15 mL of nuclease-free water.
- 10 Prepare an Eppendorf twin.tec source plate labeled *Wash #1*. For each well to be processed, add 160 μ L of SureSelect Wash Buffer #1.
- 11 Prepare a Nunc DeepWell source plate labeled *Wash #2*. For each well to be processed, add 800 μ L of SureSelect Wash Buffer #2.
- 12 Place the silver Nunc DeepWell plate insert on position 6 of the Bravo deck. This insert is required to facilitate heat transfer to DeepWell source plate wells during the Capture&Wash runset.
- 13 Place the *Wash #2* source plate on the silver insert at position 6 of the Bravo deck. Make sure the plate is seated properly on the silver DeepWell insert.

Load the Agilent NGS Workstation

14 Load the Labware MiniHub according to [Table 58](#), using the plate orientations shown in [Figure 4](#).

Table 58 Initial MiniHub configuration for SureSelectCapture&Wash_v1.0.rst

Vertical Shelf Position	Cassette 1	Cassette 2	Cassette 3	Cassette 4
Shelf 5 (Top)	Empty	Empty	Empty	Empty
Shelf 4	Empty	Empty	Empty	Empty
Shelf 3	Empty Eppendorf plate	Empty	Wash #1 Eppendorf source plate	Empty
Shelf 2	Empty	Nuclease-free water reservoir	Empty	Empty
Shelf 1 (Bottom)	Empty	Empty	Empty	Empty tip box

15 Load the Bravo deck according to [Table 59](#) (positions 5 and 6 should already be loaded).

Table 59 Initial Bravo deck configuration for SureSelectCapture&Wash_v1.0.rst

Location	Content
1	Empty waste reservoir (Axygen 96 Deep Well Plate, square wells)
4	Empty red aluminum insert
5	Dynabeads M-270 streptavidin bead DeepWell source plate
6	Wash #2 DeepWell source plate seated on silver insert

4 Hybridization

Step 3. Capture the hybridized DNA

16 Load the BenchCel Microplate Handling Workstation according to Table 60.

Table 60 Initial BenchCel configuration for SureSelectCapture&Wash_v1.0.rst

No. of Columns Processed	Rack 1	Rack 2	Rack 3	Rack 4
1	1 Tip box	Empty	Empty	Empty
2	2 Tip boxes	Empty	Empty	Empty
3	3 Tip boxes	Empty	Empty	Empty
4	4 Tip boxes	Empty	Empty	Empty
6	6 Tip boxes	Empty	Empty	Empty
12	11 Tip boxes	Empty	Empty	Empty

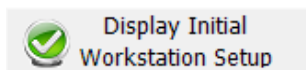
Run VWorks runset SureSelectCapture&Wash_v1.0.rst

17 On the SureSelect setup form, under **Select Protocol to Run**, select **SureSelectCapture&Wash_v1.0.rst**.

18 Under **Select PCR plate labware for Thermal Cycling**, select the specific type of PCR plate containing the DNA samples at position 6.

19 Select the number of columns of samples to be processed. Runs must include 1, 2, 3, 4, 6, or 12 columns.

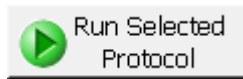
20 Click **Display Initial Workstation Setup**.



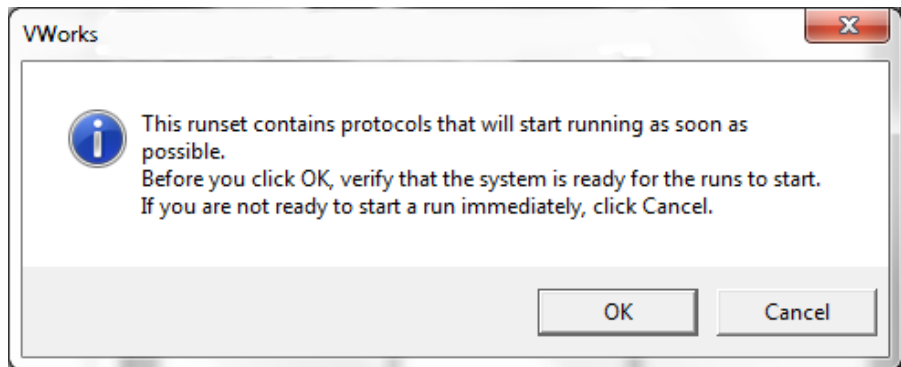
21 Verify that the NGS workstation has been set up as displayed in the Workstation Setup region of the form.



22 When verification is complete, click **Run Selected Protocol**.



23 When ready to begin the run, click **OK** in the following window. If the temperature of Bravo deck position 4 was not pre-set to 66°C, the runset will pause while position 4 reaches temperature.



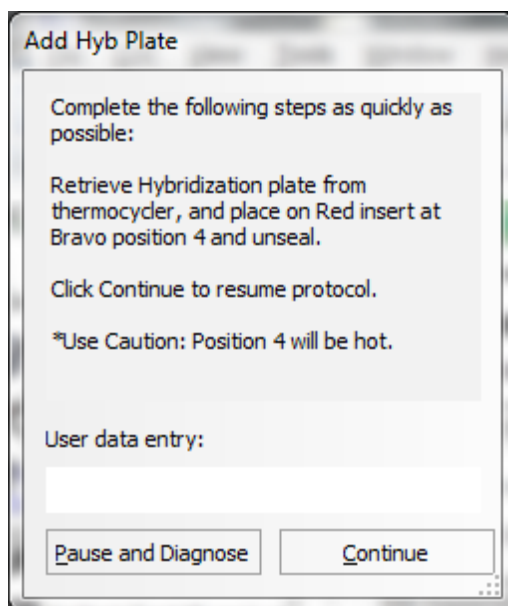
4 Hybridization

Step 3. Capture the hybridized DNA

CAUTION

It is important to complete [step 24](#) quickly and carefully. Transfer the sample plate to the Bravo platform quickly to retain the 65°C sample temperature. Unseal the plate without tilting or jerking the plate to avoid sample splashing. Make sure that the Agilent NGS Workstation is completely prepared, with deck platforms at temperature and all components in place, before you transfer the sample plate to the Bravo deck.

24 When prompted by VWorks as shown below, quickly remove the PCR plate, containing the hybridization reactions held at 65°C, from the thermal cycler. Unseal the plate carefully to avoid splashing, and quickly transfer the plate to position 4 of the Bravo deck, seated in red insert. Click **Continue** to resume the runset.

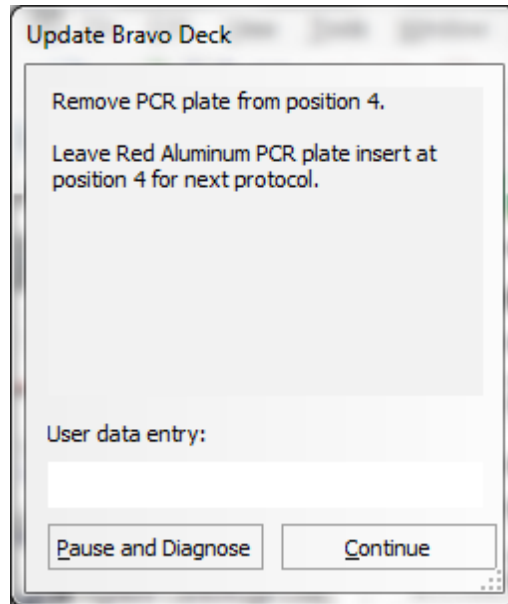


WARNING

Bravo deck position 4 will be hot.

Use caution when handling components that contact heated deck positions.

- 25** When the hybridization samples have been transferred from the PCR plate to the capture plate wells, you will be prompted by VWorks as shown below. Remove the PCR plate from position 4 of the Bravo deck, leaving the red insert in place. When finished, click **Continue** to resume the runset.



The remainder of the SureSelectCapture&Wash_v1.0.rst runset takes approximately 1.5 hours. Once the runset is complete, the captured, bead-bound DNA samples are located in the Eppendorf plate at position 9 of the Bravo deck.

When the runset is complete, seal the plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 1.0 sec and store the plate on ice while setting up the next automation protocol.

NOTE

Captured DNA is retained on the streptavidin beads during the post-capture amplification step.

4 Hybridization

Step 3. Capture the hybridized DNA



5 Indexing and Sample Prep for Multiplexed Sequencing

- Step 1. Amplify the captured libraries to add index tags 110
- Step 2. Purify the amplified indexed libraries using AMPure XP beads 119
- Step 3. Assess DNA quality and quantity 123
- Step 4. Pool samples for multiplexed sequencing 127
- Step 5. Prepare and analyze sequencing samples 129

This chapter describes the steps to add index tags by amplification, then to purify and assess quality and quantity of the libraries in order to pool indexed samples for multiplexed sequencing.



5 Indexing and Sample Prep for Multiplexed Sequencing

Step 1. Amplify the captured libraries to add index tags

Step 1. Amplify the captured libraries to add index tags

In this step, the Agilent NGS Workstation completes the liquid handling steps for PCR-based addition of indexing tags to the target-enriched DNA samples. After the PCR plate is prepared by the Agilent NGS Workstation, you transfer the plate to a thermal cycler for amplification.

CAUTION

To avoid cross-contaminating libraries, set up PCR master mixes in a dedicated clean area or PCR hood with UV sterilization and positive air flow.

Assign indexes to DNA samples

Select the appropriate indexing primer for each sample.

Use a different index primer for each sample to be sequenced in the same lane. The number of samples that may be combined per lane depends on the sequencing platform performance and the probe design size. See [Table 61](#) for sequence data requirement guidelines. Calculate the number of indexes that can be combined per lane based on these guidelines.

Table 61 Sequencing data requirement guidelines

Probe Size	Recommended Amount of Sequencing Data per Sample
1 kb up to 499 kb	0.1 to 50 Mb*
0.5 Mb up to 2.9 Mb	50 to 290 Mb*
3 Mb up to 5.9 Mb	300 to 590 Mb*

* For custom probes, Agilent recommends analyzing 100X amount of sequencing data compared to the probe design size for each sample. Pool samples according to your expected sequencing output.

Prepare the workstation

- 1 Clear the Labware MiniHub and BenchCel of plates and tip boxes.
- 2 Gently wipe down the Labware MiniHub, Bravo decks, and BenchCel with a DNA Away decontamination wipe.
- 3 Turn on the ThermoCube, set to 0°C, at position 9 of the Bravo deck. Be sure that the chiller reservoir contains at least 300 mL of 25% ethanol. Place the silver insert at position 9.
- 4 Pre-set the temperature of Bravo deck positions 4 and 6 to 4°C using the Inheco Multi TEC control touchscreen, as described in [Setting the Temperature of Bravo Deck Heat Blocks](#). On the control touchscreen, Bravo deck positions 4 corresponds to CPAC 2, position 1, while deck position 6 corresponds to CPAC 2, position 2.

5 Indexing and Sample Prep for Multiplexed Sequencing

Step 1. Amplify the captured libraries to add index tags

Prepare the index and PCR Master Mix source plates

- 5 Add 5 μL of the appropriate indexing primer to the appropriate wells of a PCR plate.

The well position for each index should correspond to the position of the RNA sample assigned to that index in the original total RNA sample plate. Keep the plate on ice.

- 6 Prepare the appropriate volume of Post-capture PCR master mix, according to [Table 62](#). Mix well using a vortex mixer and keep on ice.

CAUTION

The PCR Master Mix used at this step is highly viscous and thorough mixing prior to use is critical for optimal kit performance. Mix by vortexing at high speed for 5 seconds before removing an aliquot for use and after combining with other solutions. Pipetting up and down is not sufficient to mix this reagent.

Table 62 Preparation of Post-capture PCR Master Mix for Post-CapturePCR_RNASeq_ILM_v1.0.pro

SureSelect ^{XT} Reagent	Volume for 1 Library	Volume for 1 Column	Volume for 2 Columns	Volume for 3 Columns	Volume for 4 Columns	Volume for 6 Columns	Volume for 12 Columns
PCR Master Mix* (bottle)	25.0 μL	307.5 μL	512.5 μL	717.5 μL	922.5 μL	1332.5 μL	2665 μL
ILM Post-Capture PCR Primer† (green cap)	1.0 μL	12.3 μL	20.5 μL	28.7 μL	36.9 μL	53.3 μL	106.6 μL
Total Volume	26.0 μL	319.8 μL	533.0 μL	746.2 μL	959.4 μL	1385.8 μL	2771.6 μL

* May also be labeled as *RNA Seq PCR Master Mix*.

† May also be labeled as *RNA Seq ILM Post-Capture PCR Primer*.

- 7 Using the same Nunc DeepWell master mix source plate that was used for the SureSelectHybridization_v1.0.pro protocol, add the volume of Post-capture PCR master mix indicated in [Table 63](#) to all wells of column 4 of the master mix source plate. The final configuration of the master mix source plate is shown in [Figure 12](#).

Table 63 Preparation of the Master Mix Source Plate for Post-CapturePCR_RNASeq_ILM_v1.0.pro

Master Mix Solution	Position on Source Plate	Volume of Master Mix added per Well of Nunc Deep Well Source Plate					
		1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
Post-capture PCR Master Mix (from Table 62)	Column 4 (A4-H4)	36.7 µL	63.4 µL	90.0 µL	116.7 µL	170.0 µL	343.2 µL

NOTE

If you are using a new DeepWell plate for the post-capture PCR source plate (for example, when amplifying the second half of the captured DNA sample), leave columns 1 to 3 empty and add the PCR Master Mix to column 4 of the new plate.

5 Indexing and Sample Prep for Multiplexed Sequencing

Step 1. Amplify the captured libraries to add index tags

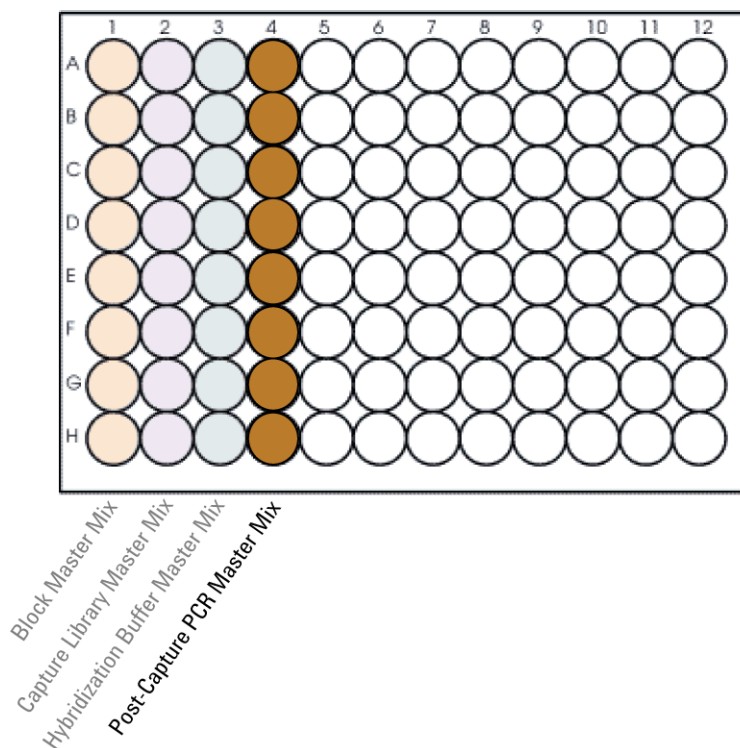


Figure 12 Configuration of the master mix source plate for Post-CapturePCR_RNaseq_ILM_v1.0.pro. Columns 1–3 were used to dispense master mixes for the SureSelectHybridization_v1.0.pro protocol.

- 8 Seal the master mix source plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 1.0 sec.
- 9 Centrifuge the plate for 30 seconds to drive the well contents off the walls and plate seal and to eliminate any bubbles.

Load the Agilent NGS Workstation

10 Load the Labware MiniHub according to [Table 64](#), using the plate orientations shown in [Figure 4](#).

Table 64 Initial MiniHub configuration for Post-CapturePCR_RNASeq_ILM_v1.0.pro

Vertical Shelf Position	Cassette 1	Cassette 2	Cassette 3	Cassette 4
Shelf 5 (Top)	Empty	Empty	Empty	Empty
Shelf 4	Empty	Empty	Empty	Empty
Shelf 3	Empty	Empty	Empty	Empty
Shelf 2	Empty tip box	Empty	Empty	Empty
Shelf 1 (Bottom)	New tip box	Empty	Empty	Empty tip box

11 Load the Bravo deck according to [Table 65](#).

Table 65 Initial Bravo deck configuration for Post-CapturePCR_RNASeq_ILM_v1.0.pro

Location	Content
4	Captured DNA bead suspensions in Eppendorf twin.tec plate
6	Indexing primers in PCR plate seated on red insert (PCR plate type must be specified on setup form under step 2)
9	Master mix plate containing Post-capture PCR Master Mix in Column 4 (Nunc DeepWell plate seated on silver insert)

5 Indexing and Sample Prep for Multiplexed Sequencing

Step 1. Amplify the captured libraries to add index tags

12 Load the BenchCel Microplate Handling Workstation according to Table 66.

Table 66 Initial BenchCel configuration for Post-CapturePCR_RNASeq_ILM_v1.0.pro

No. of Columns Processed	Rack 1	Rack 2	Rack 3	Rack 4
1	1 Tip box	Empty	Empty	Empty
2	1 Tip box	Empty	Empty	Empty
3	1 Tip box	Empty	Empty	Empty
4	1 Tip box	Empty	Empty	Empty
6	1 Tip box	Empty	Empty	Empty
12	1 Tip box	Empty	Empty	Empty

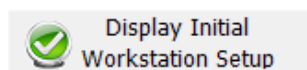
Run VWorks protocol Post-CapturePCR_RNASeq_ILM_v1.0.pro

13 On the SureSelect setup form, under **Select Protocol to Run**, select **Post-CapturePCR_RNASeq_ILM_v1.0.pro**.

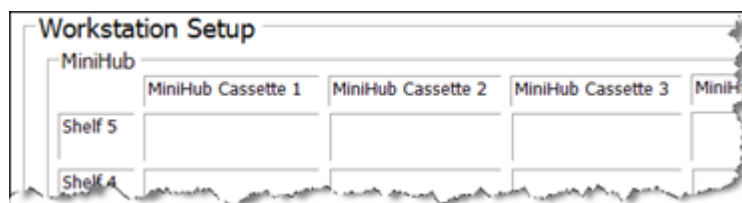
14 Under **Select PCR plate labware for Thermal Cycling**, select the specific type of PCR plate containing the indexing primers at position 6.

15 Select the number of columns of samples to be processed. Runs must include 1, 2, 3, 4, 6, or 12 columns.

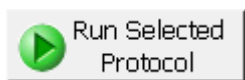
16 Click **Display Initial Workstation Setup**.



17 Verify that the NGS workstation has been set up as displayed in the Workstation Setup region of the form.

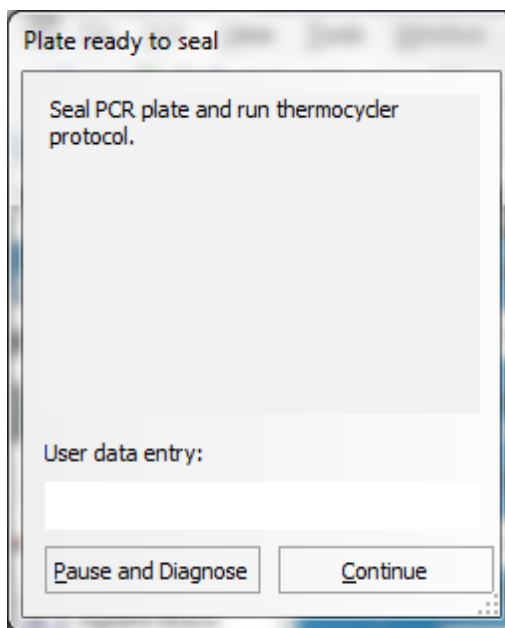


18 When verification is complete, click **Run Selected Protocol**.



Running the Post-CapturePCR_RNASeq_ILM_v1.0.pro protocol takes approximately 15 minutes. Once complete, the PCR-ready samples, containing captured DNA, indexing primer, and PCR master mix are located in the PCR plate at position 6 of the Bravo deck. The Eppendorf plate containing the remaining bead-bound captured DNA samples, which may be stored for future use at 4°C overnight, or at -20°C for longer-term storage, is located at position 4 of the Bravo deck.

19 When you see the following prompt, remove the PCR plate from position 6 of the Bravo deck and seal the plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 3.0 seconds.



5 Indexing and Sample Prep for Multiplexed Sequencing

Step 1. Amplify the captured libraries to add index tags

20 Transfer the PCR plate to a thermal cycler and run the PCR amplification program shown in [Table 67](#).

Table 67 Post-Capture PCR cycling program

Segment	Number of Cycles	Temperature	Time
1	1	95°C	2 minutes
2	12	95°C	30 seconds
		57°C	30 seconds
		72°C	1 minute
3	1	72°C	5 minutes
4	1	4°C	Hold

Step 2. Purify the amplified indexed libraries using AMPure XP beads

In this step, the Agilent NGS Workstation transfers AMPure XP beads to the indexed DNA sample plate and then collects and washes the bead-bound DNA.

Prepare the workstation and reagents

- 1 Clear the Labware MiniHub and BenchCel of all plates and tip boxes.
- 2 Gently wipe down the Labware MiniHub, Bravo decks, and BenchCel with a DNA Away decontamination wipe.
- 3 Let the AMPure XP beads come to room temperature for at least 30 minutes. *Do not freeze the beads at any time.*
- 4 Mix the bead suspension well so that the reagent appears homogeneous and consistent in color.
- 5 Prepare a Nunc DeepWell source plate containing AMPure XP beads. For each well to be processed, add 95 μL of homogeneous AMPure XP beads per well to the Nunc DeepWell plate.
- 6 Prepare a Thermo Scientific reservoir containing 15 mL of nuclease-free water.
- 7 Prepare a separate Thermo Scientific reservoir containing 45 mL of freshly-prepared 70% ethanol.

5 Indexing and Sample Prep for Multiplexed Sequencing

Step 2. Purify the amplified indexed libraries using AMPure XP beads

- 8 Load the Labware MiniHub according to [Table 68](#), using the plate orientations shown in [Figure 4](#).

Table 68 Initial MiniHub configuration for AMPureXP_v1.1.pro:Post-Capture PCR

Vertical Shelf Position	Cassette 1	Cassette 2	Cassette 3	Cassette 4
Shelf 5 (Top)	Empty Nunc DeepWell plate	Empty	Empty	Empty
Shelf 4	Empty	Empty	Empty	Empty
Shelf 3	Empty	Empty Eppendorf twin.tec plate	Empty	Empty
Shelf 2	Empty	Nuclease-free water reservoir from step 6	AMPure XP beads in Nunc DeepWell plate from step 5	Empty
Shelf 1 (Bottom)	Empty	70% ethanol reservoir from step 7	Empty	Empty tip box

- 9 Load the Bravo deck according to [Table 69](#).

Table 69 Initial Bravo deck configuration for AMPureXP_v1.1.pro:Post-Capture PCR

Location	Content
1	Empty waste reservoir (Axygen 96 Deep Well Plate, square wells)
9	Indexed library samples in PCR plate seated on red insert (PCR plate type must be specified on setup form under step 2)

Step 2. Purify the amplified indexed libraries using AMPure XP beads

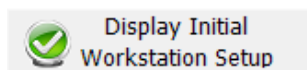
10 Load the BenchCel Microplate Handling Workstation according to Table 70.

Table 70 Initial BenchCel configuration for AMPureXP_v1.1.pro:Post-Capture PCR

No. of Columns Processed	Rack 1	Rack 2	Rack 3	Rack 4
1	1 Tip box	Empty	Empty	Empty
2	1 Tip box	Empty	Empty	Empty
3	2 Tip boxes	Empty	Empty	Empty
4	2 Tip boxes	Empty	Empty	Empty
6	3 Tip boxes	Empty	Empty	Empty
12	6 Tip boxes	Empty	Empty	Empty

Run VWorks protocol AMPureXP_v1.1.pro:Post-Capture PCR

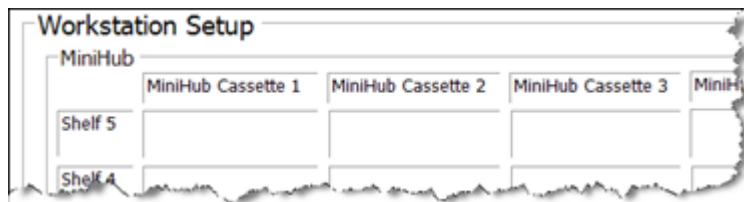
- 11** On the SureSelect setup form, under **Select Protocol to Run**, select **AMPureXP_v1.1.pro:Post-Capture PCR**.
- 12** Under **Select PCR plate labware for Thermal Cycling**, select the specific type of PCR plate containing the DNA samples at position 9.
- 13** Select the number of columns of samples to be processed. Runs must include 1, 2, 3, 4, 6, or 12 columns.
- 14** Click **Display Initial Workstation Setup**.



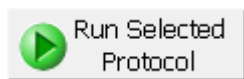
5 Indexing and Sample Prep for Multiplexed Sequencing

Step 2. Purify the amplified indexed libraries using AMPure XP beads

- 15 Verify that the NGS workstation has been set up as displayed in the Workstation Setup region of the form.



- 16 When verification is complete, click **Run Selected Protocol**.



The purification protocol takes approximately 45 minutes. When complete, the purified DNA samples are in the Eppendorf plate located on Bravo deck position 7.

Step 3. Assess DNA quality and quantity

Option 1: Analysis using the Agilent 2100 Bioanalyzer and High Sensitivity DNA Assay

See the *High Sensitivity DNA Kit Guide* at www.genomics.agilent.com for more information on doing this step.

- 1 Set up the 2100 Bioanalyzer as instructed in the High Sensitivity DNA Assay kit guide.

NOTE

Version B.02.07 or higher of the Agilent 2100 Expert Software is required for High Sensitivity DNA Assay Kit runs.

- 2 Seal the sample plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 1.0 sec.
- 3 Vortex the plate to mix samples in each well, then centrifuge the plate for 30 seconds to drive the well contents off the walls and plate seal.
- 4 Prepare the chip, samples and ladder as instructed in the reagent kit guide, using 1 µL of each sample for the analysis.

NOTE

For some samples, Bioanalyzer results are improved by diluting 1 µL of the sample in 9 µL of 10 mM Tris, 1 mM EDTA prior to analysis. Be sure to mix well by vortexing at 2000 rpm on the IKA vortex supplied with the Bioanalyzer before analyzing the diluted samples.

- 5 Load the prepared chip into the 2100 Bioanalyzer and start the run within five minutes after preparation.

Verify that the electropherogram shows an average DNA fragment size of approximately 200 to 700 bp. A sample electropherogram is shown in [Figure 13](#).

Stopping Point If you do not continue to the next step, seal the plate and store at -20°C.

5 Indexing and Sample Prep for Multiplexed Sequencing

Step 3. Assess DNA quality and quantity

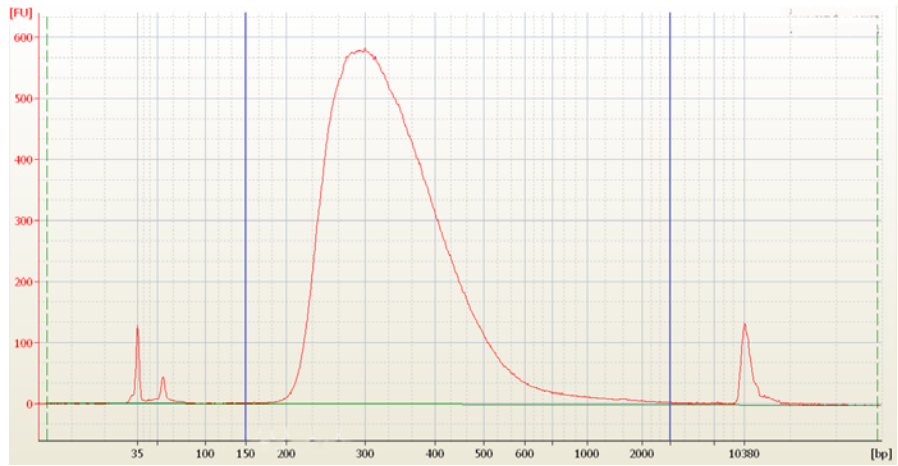


Figure 13 Analysis of indexed DNA using the High Sensitivity DNA Assay.

Option 2: Analysis using the Agilent 4200 TapeStation and High Sensitivity D1000 ScreenTape

Use a High Sensitivity D1000 ScreenTape and reagent kit to analyze the indexed DNA. For more information to do this step, see the [Agilent High Sensitivity D1000 Assay Quick Guide](#).

- 1 Seal the DNA sample plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 1.0 sec.
- 2 Vortex the plate to mix samples in each well, then centrifuge the plate for 30 seconds to drive the well contents off the walls and plate seal.
- 3 Prepare the TapeStation samples as instructed in the instrument user manual. Use 2 µL of each indexed DNA sample diluted with 2 µL of High Sensitivity D1000 sample buffer for the analysis.

CAUTION

For accurate quantitation, make sure to thoroughly mix the combined DNA and sample buffer by vortexing the assay plate or tube strip for 1 minute on the IKA MS3 vortex mixer provided with the 4200 TapeStation system before loading the samples.

If no IKA MS3 vortex mixer is available, ensure that the samples are mixed thoroughly by vortexing for 10 seconds at high speed on a manual vortex mixer.

- 4 Load the sample plate or tube strips from [step 3](#), the High Sensitivity D1000 ScreenTape, and loading tips into the TapeStation as instructed in the reagent kit guide. Start the run.
- 5 For each sample, measure the concentration of the library (ng/µL) by integrating under the peak at approximately 200 to 700 bp. A sample electropherogram is shown in [Figure 14](#).

Stopping Point If you do not continue to the next step, seal the plate and store at -20°C.

5 Indexing and Sample Prep for Multiplexed Sequencing

Step 3. Assess DNA quality and quantity

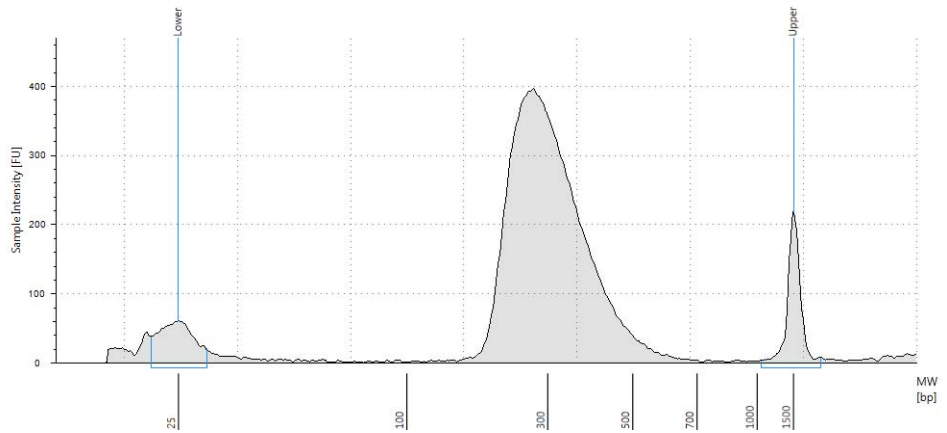


Figure 14 Analysis of purified indexed DNA amplicons using the High Sensitivity D1000 ScreenTape.

Step 4. Pool samples for multiplexed sequencing

The number of indexed libraries that may be multiplexed in a single sequencing lane is determined by the output specifications of the platform used, together with the amount of sequencing data required for your research design. Calculate the number of indexes that can be combined per lane, according to the capacity of your platform and the amount of sequencing data required per sample.

- 1 Combine the libraries such that each index-tagged sample is present in equimolar amounts in the pool. For each library, use the formula below to determine the amount of indexed sample to use.

$$\text{Volume of Index} = \frac{V(f) \times C(f)}{\# \times C(i)}$$

where $V(f)$ is the final desired volume of the pool,

$C(f)$ is the desired final concentration of all the DNA in the pool

$\#$ is the number of indexes, and

$C(i)$ is the initial concentration of each indexed sample.

Table 71 shows an example of the amount of 4 index-tagged samples (of different concentrations) and Low TE needed for a final volume of 20 μL at 10 nM.

Table 71 Example of indexed sample volume calculation for total volume of 20 μL

Component	V(f)	C(i)	C(f)	#	Volume to use (μL)
Sample 1	20 μL	20 nM	10 nM	4	2.5
Sample 2	20 μL	10 nM	10 nM	4	5
Sample 3	20 μL	17 nM	10 nM	4	2.9
Sample 4	20 μL	25 nM	10 nM	4	2
Low TE					7.6

- 2 Adjust the final volume of the pooled library to the desired final concentration.

5 Indexing and Sample Prep for Multiplexed Sequencing

Step 4. Pool samples for multiplexed sequencing

- If the final volume of the combined index-tagged samples is less than the desired final volume, $V(f)$, add Low TE to bring the volume to the desired level.
 - If the final volume of the combined index-tagged samples is greater than the final desired volume, $V(f)$, lyophilize and reconstitute to the desired volume.
- 3** If you store the library before sequencing, add Tween 20 to 0.1% v/v and store at -20°C short term.

Exact library pool dilution and processing can vary based on the flow cell capacity and analysis pipeline versions being used. Refer to the appropriate Illumina user guide for instructions.

Step 5. Prepare and analyze sequencing samples

Proceed to cluster amplification using the appropriate Illumina Paired-End Cluster Generation Kit. See [Table 72](#) for kit configurations compatible with the recommended read length.

The optimal seeding concentration for SureSelect^{XT} target-enriched RNA sequencing libraries varies according to sequencing platform, run type, and Illumina kit version. See [Table 72](#) for guidelines. Seeding concentration and cluster density may also need to be optimized based on the fragment size range for the library and on the desired output and data quality.

Follow Illumina's recommendation for a PhiX control in a low-concentration spike-in for improved sequencing quality control.

Table 72 Illumina Kit Configuration Selection Guidelines

Platform	Run Type	Read Length	SBS Kit Configuration	Chemistry	Seeding Concentration
HiSeq 2500	Rapid Run	2 × 100 bp	200 Cycle Kit	v2	10–13 pM
HiSeq 2500	High Output	2 × 100 bp	200 Cycle Kit	v3	10–13 pM
HiSeq 2500	High Output	2 × 100 bp	250 Cycle Kit	v4	13–16 pM
HiSeq 2000	All Runs	2 × 100 bp	200 Cycle Kit	v3	7–11 pM
HiSeq 2000	All Runs	2 × 100 bp	250 Cycle Kit	v4	10–14 pM
MiSeq	All Runs	2 × 100 bp	300 Cycle Kit	v2	10–13 pM
MiSeq	All Runs	2 × 75 bp	150 Cycle Kit	v3	14–19 pM
NextSeq 500/550	All Runs	2 × 100 bp	300 Cycle Kit	v2	1.7–2.0 pM
HiSeq 3000/4000	All Runs	2 × 100 bp	300 Cycle Kit	v1	200 pM

5 Indexing and Sample Prep for Multiplexed Sequencing

Step 5. Prepare and analyze sequencing samples

Sequencing run setup guidelines for 8-bp indexes

Sequencing runs must be set up to perform an 8-bp index read. See the [Reference](#) chapter for complete index sequence information.

For the HiSeq and NextSeq 500 (v1) platforms, use the *Cycles* settings shown in [Table 73](#). Cycle number settings can be specified on the *Run Configuration* screen of the instrument control software interface after choosing *Custom* from the index type selection buttons.

Table 73 Cycle Number settings for HiSeq and NextSeq platforms

Run Segment	Cycle Number
Read 1	100
Index 1 (i7)	8
Index 2 (i5)	0
Read 2	100

For the MiSeq platform, use the Illumina Experiment Manager (IEM) software to generate a Sample Sheet that includes the run parameters specified in [Table 74](#).

Table 74 Run parameters for MiSeq platform Sample Sheet

Parameter	Entry
Workflow	GenerateFASTQ
Cycles for Read 1	100 for v2 chemistry 75 for v3 chemistry
Cycles for Read 2	100 for v2 chemistry 75 for v3 chemistry
Index 1 (i7) Sequence (enter in Data Section for each sample)	Type the 8-nt index sequence for each individual sample (see Table 37 on page 67).

Sequence analysis guidelines

The SureSelect^{XT} RNA sequencing library preparation method preserves RNA strandedness using dUTP second-strand marking. The sequence of read 1, which starts at the P5 end, matches the reverse complement of the poly-A RNA transcript strand. Read 2, which starts at the P7 end, matches the poly-A RNA transcript strand. When running analysis of this data to determine strandedness, it is important to include this information. For example, when using the Picard tools (<https://broadinstitute.github.io/picard>) to calculate RNA sequencing metrics, it is important to include the parameter *STRAND_SPECIFICITY=SECOND_READ_TRANSCRIPTION_STRAND* to correctly calculate the strand specificity metrics.

5 Indexing and Sample Prep for Multiplexed Sequencing

Step 5. Prepare and analyze sequencing samples



6 Reference

Kit Contents 134

Nucleotide Sequences of SureSelect^{XT} Indexes A01 to H12 138

Plasticware quantities for automation protocols 139

This chapter contains reference information, including component kit contents, index sequences, and plasticware requirements.



Kit Contents

The SureSelect^{XT} RNA Reagent Kit includes the component kits listed in [Table 75](#). The contents of each component kit are detailed in [Table 76](#) through [Table 79](#).

Table 75 SureSelect^{XT} RNA Reagent Kit Content

Component Kits	Storage Condition	Part Number
SureSelect RNA Library Prep, ILM (Pre PCR) [*]	-20°C	5500-0135
SureSelect Poly-A Selection Module (Pre PCR) [†]	4°C	5190-6411
SureSelect Target Enrichment Box 1	Room Temperature	5190-8646
SureSelect Target Enrichment Box 2	-20°C	5190-6262

* May also be labeled as *SureSelect Strand Specific RNA Library Prep, ILM, Box 1*.

† May also be labeled as *SureSelect Strand Specific RNA Library Prep, ILM, Box 2*.

Table 76 SureSelect RNA Library Prep, ILM (Pre PCR) Content

Equivalent RNA Library Preparation Kit Component Names*		Format (96 Reaction Kit, p/n 5500-0135)
Current name	Retired name	
Fragmentation Mix	RNA Seq Fragmentation Mix	bottle
First Strand Master Mix	RNA Seq First Strand Master Mix	tube with orange cap
Second Strand Enzyme Mix [†]	RNA Seq Second-Strand + End Repair Enzyme Mix	bottle
Second Strand Oligo Mix	RNA Seq Second-Strand + End Repair Oligo Mix	tube with yellow cap
dA Tailing Master Mix	RNA Seq dA Tailing Master Mix	bottle
SureSelect Ligation Master Mix	(no change)	tube with purple cap
SureSelect Oligo Adaptor Mix	(no change)	tube with blue cap
PCR Master Mix	RNA Seq PCR Master Mix	bottle
Uracil DNA Glycosylase (UDG)	(no change)	tube with yellow cap
SureSelect Primer	(no change)	tube with brown cap
ILM Reverse PCR Primer	RNA Seq ILM Reverse PCR Primer	tube with black cap
ILM Post-Capture PCR Primer	RNA Seq ILM Post-Capture PCR Primer	tube with green cap
SureSelect ^{XT} Indexes, 8 bp [‡]	(no change)	SureSelect 8 bp Indexes A01 through H12, provided in blue 96-well plate**

* Some component names were updated in June, 2020. Formulations of the reagents supplied and protocols for use of the reagents are unchanged. All components are supported through the expiration date listed on the Certificate of Analysis.

† The *Second Strand Enzyme Mix* and *Second Strand Oligo Mix* also supply the reagents used for cDNA end repair; formulations are equivalent to vials labeled as *RNA Seq Second Strand + End Repair Enzyme Mix* and *Second Strand + End Repair Oligo Mix*, respectively.

‡ See [Table 81](#) on page 138 for index sequences.

** See [Table 80](#) on page 137 for a plate map.

Table 77 SureSelect Poly-A Selection Module (Pre PCR) Content

Kit Component	Format (96 Reaction Kit, p/n 5190-6411)*
Oligo(dT) Microparticles	bottle
Bead Binding Buffer	bottle
Bead Washing Buffer	bottle
Bead Elution Buffer	bottle

* Component kit part number 5190-6411 labeled as *SureSelect Strand Specific RNA Library Prep, ILM, Box 2* also contains a vial of Nuclease Free Water and contains reagent tubes labeled as *RNA Seq Bead Binding Buffer, RNA Seq Bead Washing Buffer, and RNA Seq Bead Elution Buffer*. Formulations of the reagents supplied and protocols for use of these reagents are unchanged. Ordering information for nuclease-free water is provided in [Table 1](#) on page 12.

Table 78 SureSelect Target Enrichment Box 1 Content

Kit Component	Format (96 Reaction Kit, p/n 5190-8646)
SureSelect Hyb 1	bottle
SureSelect Hyb 2	tube with red cap
SureSelect Hyb 4	tube with black cap
SureSelect Binding Buffer	bottle
SureSelect Wash Buffer 1	bottle
SureSelect Wash Buffer 2	bottle

Table 79 SureSelect Target Enrichment Box 2 Content

Kit Component	Format (96 Reaction Kit, p/n 5190-6262)
SureSelect Hyb 3	tube with yellow cap
SureSelect Indexing Block 1	tube with green cap
SureSelect Block 2	tube with blue cap
SureSelect Indexing Block 3	tube with brown cap
SureSelect RNase Block	tube with purple cap

Table 80 Plate map for 8bp Indexes A01 through H12 provided in blue plate in Library Prep kit p/n 5500-0135

	1	2	3	4	5	6	7	8	9	10	11	12
A	A01	A02	A03	A04	A05	A06	A07	A08	A09	A10	A11	A12
B	B01	B02	B03	B04	B05	B06	B07	B08	B09	B10	B11	B12
C	C01	C02	C03	C04	C05	C06	C07	C08	C09	C10	C11	C12
D	D01	D02	D03	D04	D05	D06	D07	D08	D09	D10	D11	D12
E	E01	E02	E03	E04	E05	E06	E07	E08	E09	E10	E11	E12
F	F01	F02	F03	F04	F05	F06	F07	F08	F09	F10	F11	F12
G	G01	G02	G03	G04	G05	G06	G07	G08	G09	G10	G11	G12
H	H01	H02	H03	H04	H05	H06	H07	H08	H09	H10	H11	H12

Nucleotide Sequences of SureSelect^{XT} Indexes A01 to H12

Each index is 8 nt in length. See [page 130](#) for sequencing run setup requirements for sequencing libraries using 8-bp indexes.

Table 81 SureSelect Indexes, for indexing primers in blue 96-well plates

Index	Sequence	Index	Sequence	Index	Sequence	Index	Sequence
A01	ATGCCTAA	A04	AACTCACC	A07	ACGTATCA	A10	AATGTTGC
B01	GAATCTGA	B04	GCTAACGA	B07	GTCTGTCA	B10	TGAAGAGA
C01	AACGTGAT	C04	CAGATCTG	C07	CTAAGGTC	C10	AGATCGCA
D01	CACTTCGA	D04	ATCCTGTA	D07	CGACACAC	D10	AAGAGATC
E01	GCCAAGAC	E04	CTGTAGCC	E07	CCGTGAGA	E10	CAACCACA
F01	GACTAGTA	F04	GCTCGGTA	F07	GTGTTCTA	F10	TGGAACAA
G01	ATTGGCTC	G04	ACACGACC	G07	CAATGGAA	G10	CCTCTATC
H01	GATGAATC	H04	AGTCACTA	H07	AGCACCTC	H10	ACAGATTC
A02	AGCAGGAA	A05	AACGCTTA	A08	CAGCGTTA	A11	CCAGTTCA
B02	GAGCTGAA	B05	GGAGAACA	B08	TAGGATGA	B11	TGGCTTCA
C02	AAACATCG	C05	CATCAAGT	C08	AGTGGTCA	C11	CGACTGGA
D02	GAGTTAGC	D05	AAGGTACA	D08	ACAGCAGA	D11	CAAGACTA
E02	CGAACTTA	E05	CGCTGATC	E08	CATACCAA	E11	CCTCCTGA
F02	GATAGACA	F05	GGTGCGAA	F08	TATCAGCA	F11	TGGTGGTA
G02	AAGGACAC	G05	CCTAATCC	G08	ATAGCGAC	G11	AACAACCA
H02	GACAGTGC	H05	CTGAGCCA	H08	ACGCTCGA	H11	AATCCGTC
A03	ATCATTCC	A06	AGCCATGC	A09	CTCAATGA	A12	CAAGGAGC
B03	GCCACATA	B06	GTACGCAA	B09	TCCGTCTA	B12	TTCACGCA
C03	ACCACTGT	C06	AGTACAAG	C09	AGGCTAAC	C12	CACCTTAC
D03	CTGGCATA	D06	ACATTGGC	D09	CCATCCTC	D12	AAGACGGA
E03	ACCTCCAA	E06	ATTGAGGA	E09	AGATGTAC	E12	ACACAGAA
F03	GCGAGTAA	F06	GTCGTAGA	F09	TCTTCACA	F12	GAACAGGC
G03	ACTATGCA	G06	AGAGTCAA	G09	CCGAAGTA	G12	AACCGAGA
H03	CGGATTGC	H06	CCGACAAC	H09	CGCATACA	H12	ACAAGCTA

Plasticware quantities for automation protocols

The tables below show the quantity of each plasticware type used in each automation protocol in the workflow. Quantities listed in the tables only include unique labware that was not used in other protocols or runsets. For example, Nunc DeepWell master mix plates may be reused in multiple protocols but are counted below only where first used.

mRNA_Purification_v1.0.pro

Labware	1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
Tip boxes (filled)	2	3	4	4	6	10
Empty tip boxes (for waste tips)	2	2	2	2	2	2
Nunc DeepWell Plates	5	5	5	5	5	5
96 Eppendorf twin.tec full-skirt plates	3	3	3	3	3	3
PCR plates (compatible with thermal cycler)	4	4	4	4	4	4
Axygen square-well plate (waste)	1	1	1	1	1	1

AMPureXP_v1.1.pro:First Strand

Labware	1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
Tip boxes (filled)	1	1	2	2	3	6
Empty tip boxes (for waste tips)	1	1	1	1	1	1
Nunc DeepWell Plates	2	2	2	2	2	2
96 Eppendorf twin.tec full-skirt plates	1	1	1	1	1	1
Thermo Scientific Reservoirs	2	2	2	2	2	2
Axygen square-well plate (waste)	1	1	1	1	1	1

6 Reference

Plasticware quantities for automation protocols

LibraryPrep_RNASeq_ILM_v1.1.rst

Labware	1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
Tip boxes (filled)	2	3	4	5	7	14
Empty tip boxes (for waste tips)	1	1	1	1	1	1
Nunc DeepWell Plates	3	3	3	3	3	3
96 Eppendorf twin.tec full-skirt plates	4	4	4	4	4	4
Thermo Scientific Reservoirs	2	2	2	2	2	2
Axygen square-well plate (waste)	1	1	1	1	1	1

Pre-CapturePCR_RNASeq_ILM_v1.0.pro

Labware	1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
Tip boxes (filled)	1	1	1	1	1	1
Empty tip boxes (for waste tips)	1	1	1	1	1	1
PCR plates (compatible with thermal cycler)	1	1	1	1	1	1

AMPureXP_v1.1.pro:Pre-Capture PCR

Labware	1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
Tip boxes (filled)	1	1	2	2	3	6
Empty tip boxes (for waste tips)	1	1	1	1	1	1
Nunc DeepWell Plates	2	2	2	2	2	2
96 Eppendorf twin.tec full-skirt plates	1	1	1	1	1	1
Thermo Scientific Reservoirs	2	2	2	2	2	2
Axygen square-well plate (waste)	1	1	1	1	1	1

Aliquot_Libraries_v1.0.pro

Labware	1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
Tip boxes (filled)	1	1	1	1	1	1
Empty tip boxes (for waste tips)	1	1	1	1	1	1
PCR plates (compatible with thermal cycler)	1	1	1	1	1	1

SureSelectHybridization_v1.0.pro

Labware	1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
Tip boxes (filled)	1	1	2	2	3	5
Empty tip boxes (for waste tips)	1	1	1	1	1	1
Nunc DeepWell Plates	1	1	1	1	1	1
96 Eppendorf twin.tec full-skirt plates	1	1	1	1	1	1

SureSelectCapture&Wash_v1.0.rst

Labware	1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
Tip boxes (filled)	1	2	3	4	6	11
Empty tip boxes (for waste tips)	1	1	1	1	1	1
Nunc DeepWell Plates	2	2	2	2	2	2
96 Eppendorf twin.tec full-skirt plates	2	2	2	2	2	2
Thermo Scientific Reservoir	1	1	1	1	1	1
Axygen square-well plate (waste)	1	1	1	1	1	1

6 Reference

Plasticware quantities for automation protocols

Post-CapturePCR_RNASEq_ILM_v1.0.pro

Labware	1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
Tip boxes (filled)	2	2	2	2	2	2
Empty tip boxes (for waste tips)	2	2	2	2	2	2
PCR plates (compatible with thermal cycler)	1	1	1	1	1	1

AMPureXP_v1.1.pro:Post-Capture PCR

Labware	1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
Tip boxes (filled)	1	1	2	2	3	6
Empty tip boxes (for waste tips)	1	1	1	1	1	1
Nunc DeepWell Plates	2	2	2	2	2	2
96 Eppendorf twin.tec full-skirt plates	1	1	1	1	1	1
Thermo Scientific Reservoirs	2	2	2	2	2	2
Axygen square-well plate (waste)	1	1	1	1	1	1

www.agilent.com

In This Book

This guide contains information to run the SureSelect^{XT} Automated RNA Target Enrichment protocol using the Agilent NGS Workstation.

© Agilent Technologies, Inc. 2013, 2014,
2017, 2018, 2020

Version G0, June 2020



G9691-90020



Agilent Technologies