

# **HaloPlex Exome Target Enrichment System**

## **Automation Protocol**

### **For Illumina Sequencing**

## **Protocol**

Version C0, July 2015



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



**Agilent Technologies**

## Notices

© Agilent Technologies, Inc. 2015

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

G9906-90020

### Edition

Version C0, July 2015

Printed in USA

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

### Technical Support

For technical product support, contact your local Agilent Support Services representative.

For US and Canada, call (800) 227-9770 (option 3,4,4). For other countries, find your support center telephone numbers at [www.agilent.com/chem/contactus](http://www.agilent.com/chem/contactus).

Or send an e-mail to:  
[SureSelect.Support@agilent.com](mailto:SureSelect.Support@agilent.com)

### Notice to Purchaser

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

## 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).

## 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 automation protocol for using the HaloPlex exome target enrichment system to prepare sequencing library samples for Illumina paired-end multiplexed sequencing platforms. Sample processing steps are automated using the Agilent NGS Bravo Option A.

### **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 Bravo for HaloPlex Exome Target Enrichment**

This chapter contains an orientation to the Agilent NGS Bravo, an overview of the HaloPlex exome target enrichment protocol, and considerations for designing HaloPlex experiments for automated processing using the Agilent NGS Bravo.

### **3 Sample Preparation**

This chapter describes the steps of the automated HaloPlex workflow to prepare target-enriched sequencing libraries for the Illumina platform.

### **4 Reference**

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

## What's New in Version C.0

- Updated product labeling statement.

## What's New in Version B.0

- Support for kits supplied with either of two indexing primer configurations.

Kits with revised index configuration (typically received December, 2014 or later) include indexing primers A01 through H12 provided in a blue plate. For kit content details see [page 92](#). For nucleotide sequences of the 8-bp indexes in this revised configuration, see [Table 39](#) on page 95.

Kits with original index configuration (typically received before December, 2014), include indexing primers 1–96 provided in a clear plate. For kit content details see [page 96](#). For nucleotide sequences of the 8-bp indexes in this original configuration, see [Table 43](#) on page 99 through [Table 48](#) on page 104.

Protocol steps for indexing using primers provided in either configuration are identical.

- Updated ordering information for Agilent 2200 TapeStation consumables ([Table 3](#) on page 11)
- Instructions for obtaining Agilent's SureCall analysis software ([page 89](#))

# Content

## 1 Before You Begin

Procedural Notes	8
Safety Notes	8
Required Reagents	9
Required Equipment	10
Optional Validation Reagents and Equipment	11
Using the Agilent NGS Workstation Option B for HaloPlex Automation	12

## 2 Using the Agilent NGS Bravo for HaloPlex Exome Target Enrichment

About the Agilent NGS Bravo Option A	14
About the Bravo Platform	14
VWorks Automation Control Software	17
Overview of the HaloPlex Exome Target Enrichment Procedure	23
Experimental Setup Considerations for Automated Runs	25
Considerations for Placement of gDNA Samples in 96-well Plates for Automated Processing	26
Considerations for Equipment Setup	26
Run Time Considerations	27

## 3 Sample Preparation

Step 1. Digest genomic DNA with restriction enzymes	30
Step 2. Hybridize HaloPlex probes A-H to DNA digests A-H for target enrichment and sample indexing	43
Step 3. Pool hybridized DNA samples	55
Step 4. Capture and amplify the target DNA	59
Step 5. Purify the amplified target DNA	73
Step 6. Remove primer dimers in second serial purification	77
Step 7. Validate enrichment and quantify enriched target DNA	81
Step 8. Pool samples with different indexes for multiplexed sequencing	85

## 4 Reference

Reference Information for Kits with Revised Index Configuration (indexing primers in blue plate) 92

Kit Contents-Revised Configuration 92

Nucleotide Sequences of HaloPlex Indexes (indexing primers in blue plate) 95

Reference Information for Kits with Original Index Configuration (indexing primers in clear plate) 96

Kit Contents-Original Configuration 96

Nucleotide Sequences of HaloPlex Indexes (indexing primers in clear plate) 99

Qualitative analysis of enrichment by gel electrophoresis 105



# 1

## Before You Begin

Procedural Notes 8

Safety Notes 8

Required Reagents 9

Required Equipment for HaloPlex Exome Target Enrichment Automated  
Protocols 10

Optional Validation Reagents and Equipment 11

Using the Agilent NGS Workstation Option B for HaloPlex Automation 12

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



## Procedural Notes

- The protocols in this manual are for use with Agilent's G9906D HaloPlex Exome Target Enrichment System for automation. The protocol is not compatible with other versions of HaloPlex reagent kits, including custom kits, Cardiomyopathy Research Panel kits, and Cancer Research Panel kits.
- The 96 reaction kit contains enough reagents to prepare master mixes for four runs of 3 columns of samples (24 samples) per run. When processing samples using runs with fewer than 24 samples, some reagents may be depleted before 96 samples are run.
- The HaloPlex protocol is optimized for digestion of 200 ng of genomic DNA (split among 8 different restriction digestion reactions) plus 25 ng excess DNA, for a total of 225 ng genomic DNA. Using lower amounts of DNA in the enrichment protocol can adversely affect your results. Use a fluorometry-based DNA quantitation method, such as PicoGreen stain or Qubit fluorometry to quantify the DNA starting material.
- Always keep pre-amplification and post-amplification DNA samples in separate work areas. Perform the enrichment procedure in the pre-amplification area. Open and store the amplified, enriched DNA samples only in the post-amplification area.
- Possible stopping points, where DNA samples may be stored between steps, are marked in the protocol. Store the samples at  $-20^{\circ}\text{C}$ , but do not subject the samples to multiple freeze/thaw cycles.
- Ensure that master mixes are thoroughly mixed, by pipetting up-and-down or by gentle vortexing, before distributing to the samples.
- 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 HaloPlex Exome Target Enrichment

Description	Vendor and part number
HaloPlex Exome Target Enrichment System Kit, ILM, 96 reactions*	Agilent, p/n G9906D
Herculase II Fusion Enzyme with dNTPs (100 mM; 25 mM for each nucleotide), 200 reactions	Agilent p/n 600677
Nuclease-free Water (not DEPC-treated)	Ambion Cat #AM9930
Agencourt AMPure XP Kit	Beckman Coulter Genomics
5 mL	p/n A63880
60 mL	p/n A63881
450 mL	p/n A63882
10 M NaOH, molecular biology grade	Sigma, p/n 72068
2 M acetic acid	Sigma, p/n A8976
10 mM Tris-HCl, pH 8.0 or 10 mM Tris-acetate, pH 8.0	General laboratory supplier
100% Ethanol, molecular biology grade	Sigma-Aldrich p/n E7023
Quant-iT dsDNA BR Assay Kit, for use with the Qubit fluorometer	
100 assays, 2-1000 ng	Life Technologies p/n Q32850
500 assays, 2-1000 ng	Life Technologies p/n Q32853

\* Kits contain enough reagents for 96 reactions total, including one or more control reactions using Enrichment Control DNA (ECD) samples. Each run of up to 96 samples should include one ECD control enrichment reaction.

## Required Equipment

**Table 2** Required Equipment for HaloPlex Exome Target Enrichment Automated Protocols

Description	Vendor and part number
Agilent NGS Bravo Option A <sup>†</sup> with VWorks software version 11.3.0.1195	Contact Agilent Automation Solutions for ordering information: Customerservice.automation@agilent.com
PlateLoc Thermal Microplate Sealer with Small Hotplate	Agilent p/n G5402#226
Robotic Pipetting Tips (Sterile, Filtered, 250 µL)	Agilent p/n 19477-022
Thermal Cycler	Agilent SureCycler 8800, p/n G8800A, 96 well plate module, p/n G8810A, and 384 well plate module, p/n G8820A, or equivalent thermal cycler <sup>‡</sup> and accessories
Eppendorf twin.tec full-skirted 96-well PCR plates	Eppendorf p/n 951020401 or 951020619
Eppendorf twin.tec full-skirted 384-well PCR plates*	Eppendorf p/n 951020702
Eppendorf twin.tec half-skirted 96-well PCR plates <sup>†</sup>	Eppendorf p/n 951020303
Thermo Scientific Reservoirs	Thermo Scientific p/n 1064156
Nunc DeepWell Plates, sterile, 1.3-mL well volume	Thermo Scientific p/n 260251
Axygen 96 Deep Well Plate, 2.2 mL, Square Well (waste reservoirs)	Axygen p/n P-2ML-SQ-C E & K Scientific p/n EK-2440
Magnetic separator ** 1.5 mL tube-compatible separator or Conical vial-compatible separator	Life Technologies DynaMag-2 magnet, p/n 12321D or equivalent DynaMag-15 magnet, p/n 12301D or equivalent
Benchtop microcentrifuge	VWR p/n 93000-196, or equivalent
Benchtop plate centrifuge	Labnet International MPS1000 Mini Plate Spinner p/n C1000, or equivalent
P10, P20, P200 and P1000 pipettes	Pipetman P10, P20, P200, P1000 or equivalent
NucleoClean Decontamination Wipes	Millipore p/n 3097
Qubit 2.0 Fluorometer	Life Technologies p/n Q32866
Qubit assay tubes	Life Technologies p/n Q32856
Vortex mixer	General laboratory supplier

\* Protocols are also compatible with Agilent NGS Workstation Option B. See [page 12](#) for more information.

† Thermal cycler must be compatible with 0.2 mL tubes.

‡ Compatible with Agilent SureCycler 8800.

\*\* Select the appropriate device based on run size. See [page 56](#) for the magnetic bead volume required for your run size.

## Optional Validation Reagents and Equipment

**Table 3** Reagents and Equipment for Optional Validation Methods

Description	Vendor and part number
2200 TapeStation Platform and Consumables	
2200 TapeStation	Agilent p/n G2964AA or G2965AA
High Sensitivity D1000 ScreenTape	Agilent p/n 5067-5584
High Sensitivity D1000 Reagents	Agilent p/n 5067-5585
2100 Bioanalyzer Platform and Consumables	
2100 Bioanalyzer Laptop Bundle	Agilent p/n G2943CA
2100 Bioanalyzer Electrophoresis Set	Agilent p/n G2947CA
High Sensitivity DNA Kit	Agilent p/n 5067-4626
Gel Electrophoresis Platform and Consumables	
XCell SureLock Mini-cell	Life Technologies p/n EI0001
Novex 6% Polyacrylamide, TBE Pre-cast Gels	Life Technologies p/n EC62655BOX
Novex TBE Running Buffer, 5X	Life Technologies p/n LC6675
Novex High-density TBE Sample Buffer, 5X	Life Technologies p/n LC6678
GelRed Nucleic Acid Stain, 3X in water	Biotium p/n 41001
DNA molecular weight markers	General laboratory supplier
UV-transilluminator	General laboratory supplier

## Using the Agilent NGS Workstation Option B for HaloPlex Automation

HaloPlex target enrichment protocols are compatible with the Agilent NGS Workstation Option B. Depending on the configuration of the system purchased, however, additional adapters may be required. Before initiating experiments, see [Table 4](#) below, and verify that the listed adapters are available for your workstation.

**Table 4** Adapter checklist for HaloPlex automation using Agilent NGS Workstation Option B

Adapter Description	Quantity Required for HaloPlex Automation	Agilent part number (single adapter)
384-well plate insert	2	G5498B#60
96-well PCR plate insert (red)*	2	G5498B#13

\* If your NGS Workstation Option B system is already equipped with one red insert, purchase one additional insert using the ordering information shown.



## 2

# Using the Agilent NGS Bravo for HaloPlex Exome Target Enrichment

About the Agilent NGS Bravo Option A 14

Overview of the HaloPlex Exome Target Enrichment Procedure 23

Experimental Setup Considerations for Automated Runs 25

This chapter contains an orientation to the Agilent NGS Bravo (Option A), an overview of the HaloPlex exome target enrichment protocol, and considerations for designing HaloPlex exome experiments for automated processing using the Agilent NGS Bravo.



## About the Agilent NGS Bravo Option A

### About the Bravo Platform

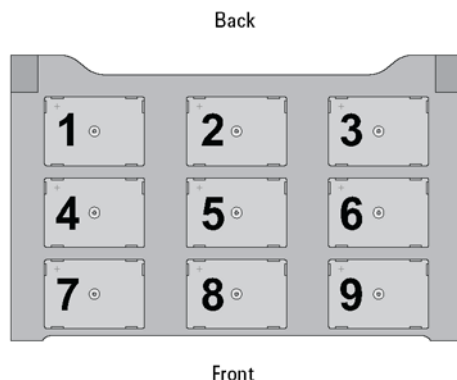
The Bravo platform is a versatile liquid handler with a nine plate-location platform deck, suitable for handling 96-well, 384-well, and 1536-well plates. The Bravo platform is controlled by the VWorks Automation Control software. Fitted with a choice of seven interchangeable fixed-tip or disposable-tip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

#### CAUTION

Before you begin, make sure that you have read and understand operating, maintenance and safety instructions for using your Bravo platform. Refer to the *Bravo Platform User Guide* (G5409-90006) and the *VWorks Software User Guide* (G5415-90063).

### 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

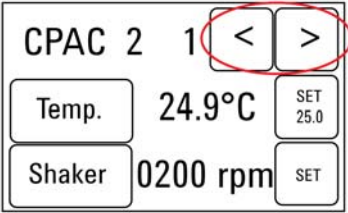
Bravo deck positions 4 and 6 are equipped with Inheco heat blocks, used to incubate sample plates at defined temperatures during the run. Runs that include low-temperature (4°C) or high-temperature (54°C) incubation steps may be expedited by pre-setting the temperature of the affected block before starting 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 5](#) for designations of the heat block-containing Bravo deck positions on the Multi TEC control device.

Table 5 Inheco Multi TEC Control touchscreen designations

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

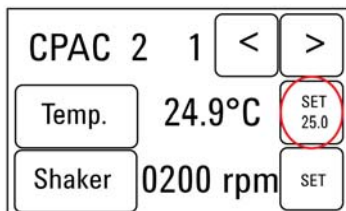
- 1 Using the arrow buttons, select the appropriate block (CPAC 2 block 1 or CPAC 2 block 2).



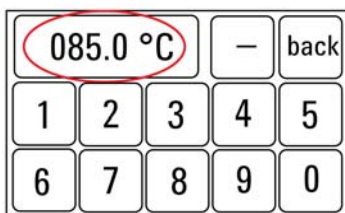
## 2 Using the Agilent NGS Bravo for HaloPlex Exome Target Enrichment

### About the Bravo Platform

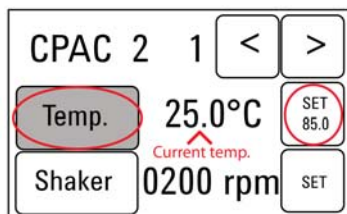
- 2 To set the temperature of the selected block, press the SET button.



- 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.





## VWorks Automation Control Software

VWorks software, included with your Agilent NGS Bravo, allows you to control the robot and integrated devices using a PC. The PC supplied with the Agilent NGS Bravo Option A is preloaded with VWorks software containing all of the necessary HaloPlex 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 HaloPlex workflow, 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 11.3.0.1195. If you have questions about VWorks version compatibility, please contact [service.automation@agilent.com](mailto:service.automation@agilent.com).

### Logging in to the VWorks software

- 1 Double-click the VWorks icon or the HaloPlex\_Exome.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.)

## 2 Using the Agilent NGS Bravo for HaloPlex Exome Target Enrichment

### VWorks Automation Control Software

#### Using the HaloPlex\_Exome.VWForm to setup and start a run

Use the VWorks form HaloPlex\_Exome.VWForm, shown below, to set up and start each HaloPlex automation protocol.

- 1 Open the form using the HaloPlex\_Exome.VWForm shortcut on your desktop.
- 2 Use the drop-down menus on the form to select the appropriate HaloPlex workflow step and number of columns of samples for the run.
- 3 Once all run parameters have been specified on the form, click **Update layout and information**.

#### NOTE

The displayed protocol will not run unless the **Update layout and information** button has been clicked.

- 4 The **Bravo Deck Setup** region of the form will then display the required placement of reaction components and labware on the Bravo deck for the specified run parameters.



## Agilent HaloPlex Automation

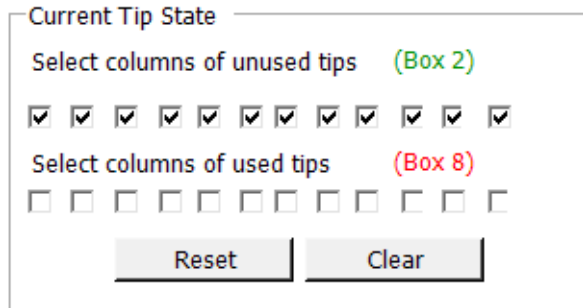


Parameters	Bravo Deck Setup	Current Tip State
1) Step: 01 Digestion_Exome.pro	New tip box	Select columns of unused tips (Box 2) [X] [X] [X] [X] [X] [X] [X] [X]
2) Number of columns of samples: 1	DNA in 96 Eppendorf Twin.tec plate on red insert 4°C	Select columns of used tips (Box 8) [X] [X] [X] [X] [X] [X] [X] [X]
3) Update layout and information	Empty 384 Eppendorf Twin.tec plate on 384 insert 4°C	Reset Clear
4) Update current tip state	Empty tip box	Reference
Status	RE Master Mixes in 96 Eppen. Twin.tec plate on red insert 0°C	Final DNA Location Labware Needs
Elapsed Time: 00:00:51		Protocol Duration Temperature Presets
Controls	Information	Advanced Settings
Start Pause Screen	It is okay to turn on the MTC to pre-chill positions 4 and 6 Turn on Thermocube to chill position 9	<input checked="" type="checkbox"/> Enable audio alerts <input type="checkbox"/> Ignore all incubation times (testing only)

C:\VWorks Workspace\NGS Option A\HaloPlex\_Exome\Protocol Files\01 Digestion\_Exomepro v2.0

- 5 Verify that the **Current Tip State** indicator on the form (shown below) matches the configuration of unused tips in the tip box at Bravo Deck position 2.

For a fresh tip box, containing 12 columns of tips, all positions of the **Current Tip State** unused tip indicator (top portion, Box 2) should be selected, as shown below. Clicking **Reset** selects all columns for position 2.



The 'Current Tip State' form contains two sections. The top section, 'Select columns of unused tips (Box 2)', has 12 checkboxes, all of which are checked. The bottom section, 'Select columns of used tips (Box 8)', has 12 checkboxes, all of which are unchecked. At the bottom of the form are two buttons: 'Reset' and 'Clear'.

Also verify that the used tip indicator (bottom portion, Box 8) matches the configuration of used tips in the tip box at Bravo Deck position 8.

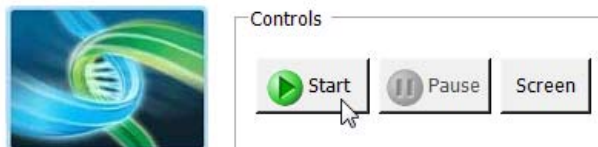
For an empty tip box, all positions of the **Current Tip State** used tip indicator (bottom portion, Box 8) should be cleared, as shown above. Clicking **Reset** clears all columns for position 8.

## NOTE

It is important that the **Current Tip State** indicator matches the configuration of tips present at Bravo Deck positions 2 and 8 when initiating the run. Tips that are inappropriately loaded onto the Bravo platform pipette head, or tips missing from the pipette head, will interfere with automated processing steps.

You can use partial tip boxes for HaloPlex automation protocols, as long as positions of available tips are accurately indicated during run setup.

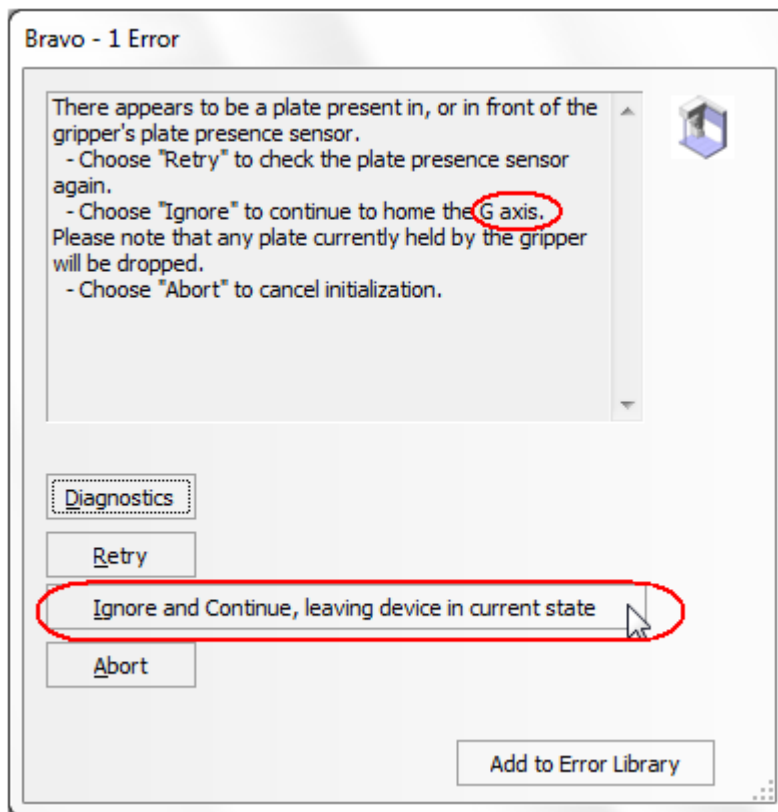
- 6 After verifying that the NGS Bravo has been set up correctly, click **Start** in the **Controls** section of the form to begin the run. Do not use the Start button on the VWorks Control Toolbar; runs must be initiated using the start button on the HaloPlex.VW Form, shown below.



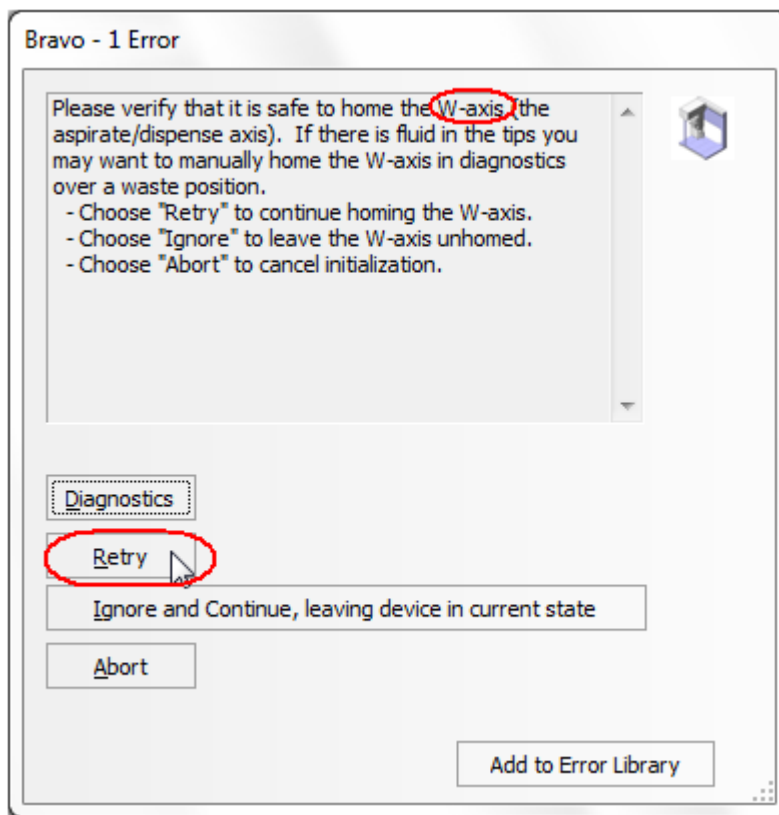
### Error messages encountered at start of run

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 Bravo 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 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 Bravo. If NGS Bravo 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 HaloPlex. VWorks form, click **Screen** in the **Controls** section of the form 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.

## Overview of the HaloPlex Exome Target Enrichment Procedure

[Figure 2](#) summarizes the HaloPlex Exome target enrichment workflow. For each sample to be sequenced, individual HaloPlex-enriched, indexed libraries are prepared. Depending on the specific sequencing platform used, multiple samples can be pooled and sequenced in a single lane.

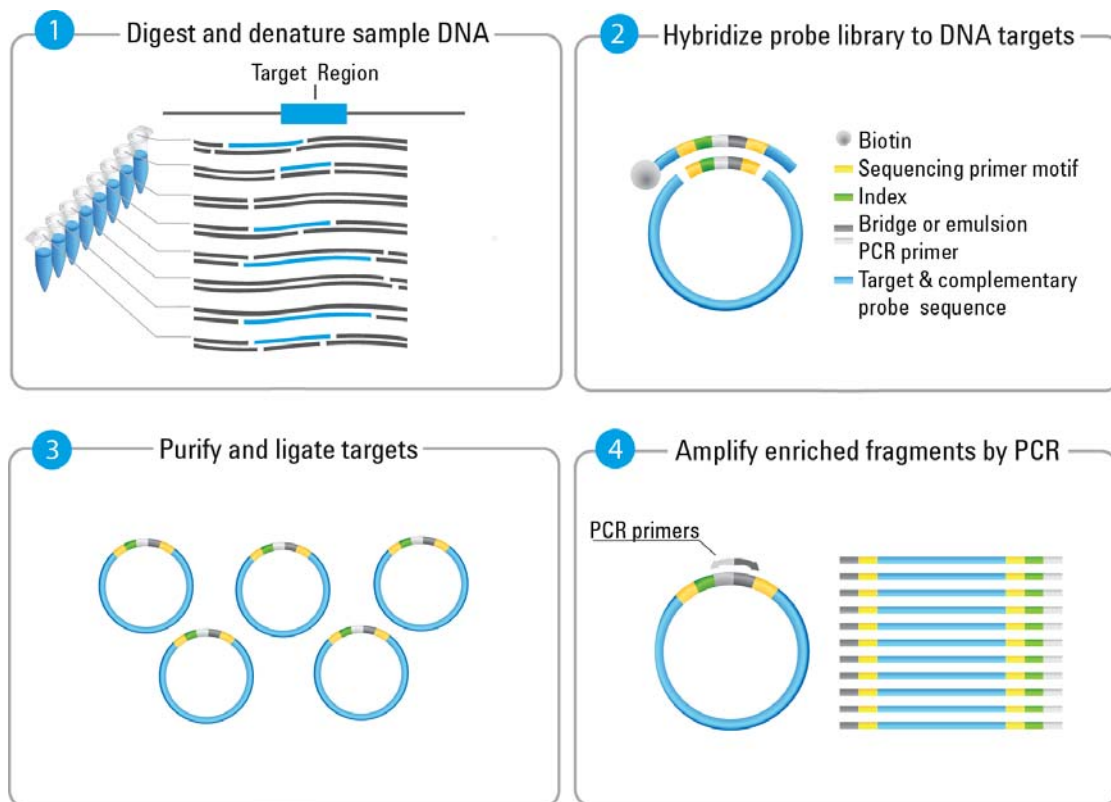
[Table 6](#) summarizes how the VWorks automation protocols are integrated into the HaloPlex Exome workflow. See the [Sample Preparation](#) chapter for complete instructions for use of the VWorks protocols for sample processing.

**Table 6** Overview of VWorks protocols used during the workflow

Workflow Step	VWorks Protocol used for Automation
Digest genomic DNA in eight separate Restriction Enzyme Master Mixes A to H	Digestion_Exome.pro
Hybridize restriction digests A to H to HaloPlex probes A to H and index DNA	Hybridization_Exome.pro
Pool hybridization reactions for each DNA sample	Post Hybridization Pooling_Exome.pro
Capture the biotin-tagged hybrids on streptavidin-coated magnetic beads and amplify the enriched DNA	Capture_Exome.pro
Purify amplified libraries	Purification_01_Exome.pro, then Purification_02_Exome.pro

## 2 Using the Agilent NGS Bravo for HaloPlex Exome Target Enrichment

### Overview of the HaloPlex Exome Target Enrichment Procedure



**Figure 2** Overall HaloPlex Exome target-enriched sequencing sample preparation workflow.



## Experimental Setup Considerations for Automated Runs

HaloPlex Automated Target Enrichment System runs may include 1, 2, 3, 4, 6, or 12 columns (equivalent to 8, 16, 24, 32, 48, or 96 wells) of gDNA samples to be enriched for sequencing on the Illumina platform. Plan your experiments using complete columns of samples.

**Table 7** 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 (see [page 92](#)) 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 gDNA Samples in 96-well Plates for Automated Processing

- The NGS Bravo processes samples column-wise beginning at column 1. gDNA 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 during hybridization to the HaloPlex probe (see [Figure 2](#)), you will need to prepare a separate plate containing the HaloPlex Indexing Primer Cassettes. Assign the wells to be indexed with their respective indexing primers during experimental design. See the [Reference](#) chapter for the nucleotide sequences of the 96 indexes used in the HaloPlex Exome Target Enrichment System.

### CAUTION

This guide includes information for kits containing two different sets of indexing primers. **Verify that you are referencing the information appropriate for your kit version before you proceed.**

Kits with indexing primers supplied in a blue plate include 8-bp indexes A01 through H12. See [page 94](#) through [page 95](#) for indexing primer A01–H12 plate map and nucleotide sequence information.

Kits with indexing primers supplied in a clear plate include 8-bp indexes 1 through 96. See [page 98](#) through [page 104](#) for indexing primer 1–96 plate map and nucleotide sequence information.

Protocol steps for indexing using primers provided in either configuration are identical.

## 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 NGS Bravo to allow rapid and efficient plate transfer.

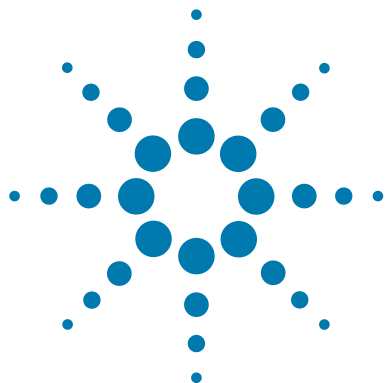
- Some workflow steps require that the sample plate be sealed, then centrifuged to collect any dispersed liquid, before being transferred between instruments. To maximize efficiency, locate the PlateLoc thermal microplate sealer and the centrifuge in close proximity to the NGS Bravo and thermal cycler.

## Run Time Considerations

The hybridization time for HaloPlex Exome enrichment is 16 to 24 hours. Based on the desired start and end time of the hybridization incubation, calculate the appropriate start time for the DNA digestion protocol, using the automation run time estimates provided in a link from the HaloPlex Exome Form in VWorks.

## **2 Using the Agilent NGS Bravo for HaloPlex Exome Target Enrichment**

### **Run Time Considerations**



### 3 Sample Preparation

- Step 1. Digest genomic DNA with restriction enzymes 30
- Step 2. Hybridize HaloPlex probes A-H to DNA digests A-H for target enrichment and sample indexing 43
- Step 3. Pool hybridized DNA samples 55
- Step 4. Capture and amplify the target DNA 59
- Step 5. Purify the amplified target DNA 73
- Step 6. Remove primer dimers in second serial purification 77
- Step 7. Validate enrichment and quantify enriched target DNA 81
- Step 8. Pool samples with different indexes for multiplexed sequencing 85

This section contains instructions for gDNA library target enrichment for sequence analysis using the Illumina platform. For each sample to be sequenced, an individual exome target-enriched, indexed library is prepared.

The HaloPlex Exome Target Enrichment System amplifies thousands of targets in the same reaction, incorporating standard Illumina paired-end sequencing motifs in the process. During hybridization, each sample can be uniquely indexed, allowing for pooling of multiple samples per sequencing lane.



## Step 1. Digest genomic DNA with restriction enzymes

In this step, gDNA samples are digested by 16 different restriction enzymes to create a library of gDNA restriction fragments. The gDNA is digested in eight different restriction reactions, each containing two restriction enzymes. **The 16 restriction enzymes are provided in two 8-well strip tubes that are distinguished by red and green color markers.** Enzymes are combined from corresponding wells of the red- and green-marked strip tubes to make eight different RE Master Mixes, which are then combined with each DNA sample in the run.

### NOTE

Successful enrichment using the protocol in this guide requires high-quality DNA samples. Before you begin, verify that the genomic DNA samples have an OD 260/280 ratio ranging from 1.8 to 2.0. Verify the size distribution of DNA in each DNA preparation by gel electrophoresis. Any smearing below 2.5 kb indicates sample degradation.

For HaloPlex target enrichment of FFPE-derived DNA samples, see Agilent publication no. G9900-90050, available at <http://www.genomics.agilent.com>. This publication provides a PCR-based protocol for assessment of DNA integrity and provides HaloPlex protocol modifications for improved performance from lower-quality DNA samples.

### Prepare the NGS Bravo

- 1 Gently wipe down the Bravo deck with a NucleoClean decontamination wipe.
- 2 Place red aluminum inserts on Bravo deck positions 4 and 9.
- 3 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.
- 4 Place a 384-well adapter insert on Bravo deck position 6.
- 5 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](#).

### NOTE

To expedite thermal cycler warm-up for the restriction digest incubation on [page 36](#), you can enter and initiate the digestion program on the thermal cycler now, and then pause the program until you are instructed to transfer the reaction plate. Be sure to pause the thermal cycler before the initiation of the 37°C incubation segment. Release the pause immediately after transferring the plate to the thermal cycler in [step 14](#) on [page 36](#). Be sure that the 384-well block is in the thermal cycler before initiating the program for warm-up.

**Prepare the DNA Sample Source Plate****NOTE**

In the protocol below, 200 ng genomic DNA is split among eight different restriction digests, with an additional 25 ng excess DNA included to allow for pipetting losses. Using <225 ng DNA in the enrichment protocol can result in low yield and can potentiate rare allele dropouts. Use a fluorometry-based DNA quantitation method, such as Qubit fluorometry or PicoGreen staining, to accurately quantify the DNA starting material.

- 1 Use the Qubit dsDNA BR Assay or PicoGreen staining kit to determine the concentration of your gDNA samples.

Follow the manufacturers instructions for the kits and instruments.

- 2 Prepare the DNA sample plate for the run, containing up to 95 gDNA samples and the Enrichment Control DNA sample, using a full-skirted 96-well Eppendorf twin.tec plate.

**NOTE**

HaloPlex Automated Target Enrichment System runs may include 1, 2, 3, 4, 6, or 12 columns of the plate. Use full columns of DNA samples for each run.

- a In well A1 of the 96-well twin.tec plate, dispense 45  $\mu$ L of the supplied Enrichment Control DNA (ECD). Store on ice.
- b In separate wells of the same 96-well twin.tec plate, dilute 225 ng of each gDNA sample in 45  $\mu$ L nuclease-free water, for a final DNA concentration of 5 ng/ $\mu$ L. Continue to store on ice.

For automated processing, fill plate wells column-wise in well order A1 to H1, then A2 to H2, ending with A12 to H12.

**Prepare the RE Master Mix Source Plate**

- 1 Prepare the appropriate amount of RE Buffer + BSA mixture, according to the table below.

**Table 8** Preparation of RE Buffer + BSA mixture for Digestion\_Exome.pro protocol

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
RE Buffer	34 $\mu$ L	408 $\mu$ L	680 $\mu$ L	952 $\mu$ L	1224 $\mu$ L	1768 $\mu$ L	3536 $\mu$ L
BSA Solution	0.85 $\mu$ L	10.2 $\mu$ L	17 $\mu$ L	23.8 $\mu$ L	30.6 $\mu$ L	44.2 $\mu$ L	88.4 $\mu$ L
<b>Total Volume</b>	<b>34.85 <math>\mu</math>L</b>	<b>418.2 <math>\mu</math>L</b>	<b>697 <math>\mu</math>L</b>	<b>975.8 <math>\mu</math>L</b>	<b>1254.6 <math>\mu</math>L</b>	<b>1812.2 <math>\mu</math>L</b>	<b>3624.4 <math>\mu</math>L</b>

### 3 Sample Preparation

#### Step 1. Digest genomic DNA with restriction enzymes

- 2 Obtain the two provided green- and red-marked Enzyme Strips from Box 1. For each strip, label the color-marked tube with A, then continue labeling the remaining tubes with B through H, in order. Keep the strips on ice.

#### CAUTION

It is important to use the restriction enzyme tube strips in the proper orientation when preparing the RE Master Mixes as described below. The red or green color marker on the tube strip and cap strip are used to mark well A of each enzyme strip.

- 3 In eight individual tubes, prepare the eight Restriction Enzyme Master Mixes A, B, C, D, E, F, G, and H according to the table below. To prepare Master Mix A, combine RE Buffer + BSA from [step 1](#) with the indicated volumes of enzyme solution from well A of the Green Enzyme Strip and from well A of the Red Enzyme Strip. Prepare Master Mixes B–H by repeating this process using enzyme solutions from the corresponding wells B–H of each provided Enzyme Strip.

**Table 9** Preparation of RE Master Mixes A–H for Digestion\_Exome.pro protocol

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
RE Buffer + BSA	4.0	51.0 µL	85.0 µL	119.0 µL	153.0 µL	221.0 µL	442.0 µL
Green Enzyme Strip enzyme A–H	0.5	6.4 µL	10.6 µL	14.9 µL	19.1 µL	27.6 µL	55.25 µL
Red Enzyme Strip enzyme A–H	0.5	6.4 µL	10.6 µL	14.9 µL	19.1 µL	27.6 µL	55.25 µL
<b>Total Volume for each Master Mix A, B, C, D, E, F, G, or H</b>	<b>5 µL</b>	<b>63.8 µL</b>	<b>106.2 µL</b>	<b>148.8 µL</b>	<b>191.2 µL</b>	<b>276.2 µL</b>	<b>552.5 µL</b>

#### NOTE

For 1-4 column runs, RE master mixes A-H may be prepared in a 8 x 0.2-mL well strip tube, using a multichannel pipette to transfer volumes from Enzyme Strips 1 and 2 to the RE master mix strip. For 6- or 12-column runs, prepare the master mixes in 1.5-mL tubes.

- 4 Mix by gentle vortexing and then spin briefly. Keep on ice.

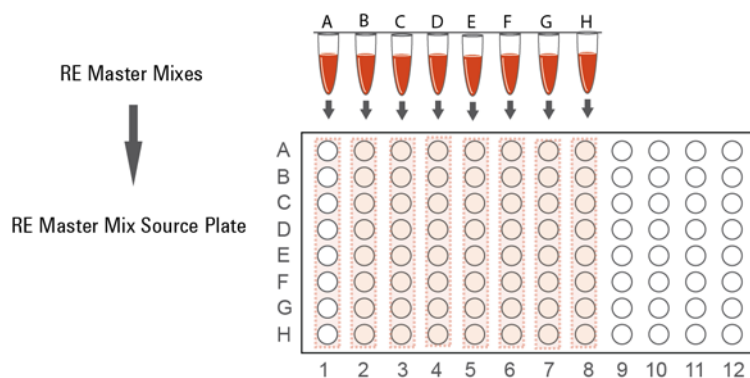


### Prepare the RE master mix source plate

- 1 Aliquot the Restriction Enzyme Master Mixes to a full-skirted 96-well Eppendorf twin.tec plate as shown in Figure 3. Add the volumes indicated in Table 10 of each master mix A–H to each well of the indicated column of the twin.tec plate. Keep the master mixes on ice during the aliquoting steps.

**Table 10** Preparation of the RE Master Mix Source Plate for Digestion\_Exome.pro

Master Mix Solution	Position on Source Plate	Volume of Master Mix added per Well of Source Plate					
		1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
RE Master Mix A	Column 1 (A1-H1)	7.3 µL	12.7 µL	18.0 µL	23.3 µL	33.9 µL	68.4 µL
RE Master Mix B	Column 2 (A2-H2)	7.3 µL	12.7 µL	18.0 µL	23.3 µL	33.9 µL	68.4 µL
RE Master Mix C	Column 3 (A3-H3)	7.3 µL	12.7 µL	18.0 µL	23.3 µL	33.9 µL	68.4 µL
RE Master Mix D	Column 4 (A4-H4)	7.3 µL	12.7 µL	18.0 µL	23.3 µL	33.9 µL	68.4 µL
RE Master Mix E	Column 5 (A5-H5)	7.3 µL	12.7 µL	18.0 µL	23.3 µL	33.9 µL	68.4 µL
RE Master Mix F	Column 6 (A6-H6)	7.3 µL	12.7 µL	18.0 µL	23.3 µL	33.9 µL	68.4 µL
RE Master Mix G	Column 7 (A7-H7)	7.3 µL	12.7 µL	18.0 µL	23.3 µL	33.9 µL	68.4 µL
RE Master Mix H	Column 8 (A8-H8)	7.3 µL	12.7 µL	18.0 µL	23.3 µL	33.9 µL	68.4 µL



**Figure 3** Preparation of the RE Master Mix source plate for automation protocol Digestion\_Exome.pro.

### 3 Sample Preparation

#### Step 1. Digest genomic DNA with restriction enzymes

##### Load the NGS Bravo and Run the Digestion\_Exome.pro VWorks Protocol

- 1 Open the HaloPlex setup form using the HaloPlex\_Exome.VWForm shortcut on your desktop.
- 2 Log in to the VWorks software.
- 3 On the setup form, under **Step**, select **01 Digestion\_Exome.pro**.

Parameters

1) Step: 01 Digestion\_Exome.pro

2) Number of columns of samples: 1

3) Update layout and information

4) Update current tip state

- 4 Select the number of columns of samples to be processed. Runs must include 1, 2, 3, 4, 6, or 12 columns.
- 5 Click **Update layout and information**.
- 6 Load the Bravo deck according to [Table 11](#).

**Table 11** Initial Bravo deck configuration for Digestion\_Exome.pro

Location	Content
1	—(empty)—
2	New tip box
3	For 12-column runs only: Empty 384-well Eppendorf twin.tec plate (no 384-well insert required) For 1- to 6-column runs: empty
4	gDNA samples in full-skirted 96-well Eppendorf twin.tec plate seated on red insert
5	—(empty)—
6	Empty 384-well Eppendorf twin.tec plate seated on 384-well insert
7	—(empty)—
8	Empty tip box
9	RE Master Mix source plate (full-skirted 96-well Eppendorf twin.tec plate) seated on red insert

- 7 Verify that the NGS Bravo has been set up as displayed in the **Bravo Deck Setup** and **Information** regions of the form.

## Step 1. Digest genomic DNA with restriction enzymes

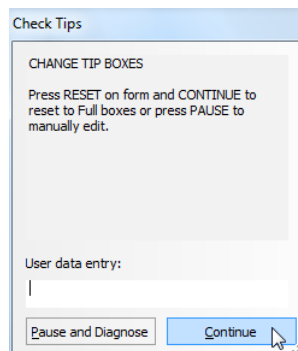
- 8 Verify that the **Current Tip State** indicator on the form matches the configuration of unused and used tips in the tip boxes at Bravo Deck positions 2 and 8, respectively. See [page 19](#) for more information on using this segment of the form during the run.
- 9 When verification is complete, click **Start** to start the run.

**NOTE**

If NGS Bravo 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 22](#) for more information.

- 10 When prompted by VWorks as shown below, replace the tip box at position 2 with a new tip box and replace the used tip box at position 8 with an empty tip box. After both tip boxes are in place, click **Reset** under **Current Tip State** on the form. Verify that the tip state was updated and then click **Continue** on the prompt shown below.

Depending on the run size, you may be prompted to change tip boxes multiple times during the run.



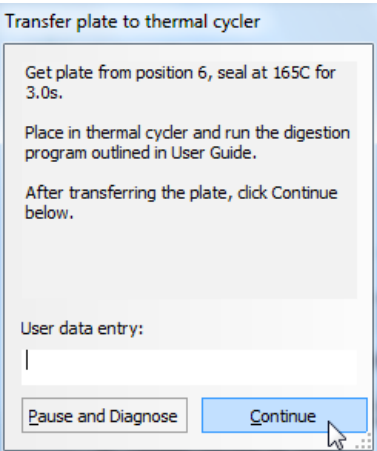
The NGS Bravo combines each gDNA sample with each RE Master Mix in wells of a 384-well reaction plate. For 1- to 6-column runs, a single 384-well restriction digest plate is prepared; for 12 column runs, two 384-well restriction digest plates are prepared.

### 3 Sample Preparation

#### Step 1. Digest genomic DNA with restriction enzymes

**11** When the NGS Bravo has finished preparing each 384-well restriction digest plate for the run, you will be prompted by VWorks as shown below.

The final Bravo deck position of the prepared restriction digest plate varies for different run sizes.



**12** Remove the 384-well plate from the Bravo deck position indicated in the prompt.

**13** Seal the sample plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 3.0 sec. Spin the plate briefly to release any bubbles trapped in the liquid.

**14** Transfer the sealed plate to a thermal cycler and run the digestion program shown in [Table 12](#), using a heated lid. After transferring the plate, click **Continue** on the prompt.

**Table 12** Thermal cycler program for HaloPlex restriction digestion

Step	Temperature	Time
Step 1	37°C	30 minutes
Step 2	8°C	Hold

Preparation of each restriction digest reaction plate takes approximately 30-45 minutes.

## Step 1. Digest genomic DNA with restriction enzymes

For 12 column-runs, the two 384-well plates are prepared sequentially, for a total run time of approximately 90 minutes. Run the thermal cycler digestion program for each plate as soon as prompted. During the 30-minute incubation of plate 1 in the thermal cycler, the NGS Bravo begins preparation of the digestion reactions in plate 2. Once the thermal cycler program is complete for plate 1, store the digested DNA in plate 1 on ice until the Digestion\_Exome.pro protocol and thermal cycler program for plate 2 is finished.

- 15** Store the 384-well DNA digestion plate(s) on ice while completing the validation steps below as well as while setting up the subsequent Hybridization protocol.

### 3 Sample Preparation

#### Step 1. Digest genomic DNA with restriction enzymes

**16** Validate the restriction digestion reaction by electrophoretic analysis of the Enrichment Control DNA (ECD) reactions.

- a** Transfer 4  $\mu$ L of each ECD digestion reaction from the wells of the 384-well reaction plate indicated in [Table 13](#) to fresh 0.2-mL PCR tubes. Note that for 12-column runs, four of the eight ECD digests are found on the first 384-well plate, and the remaining four digests are on the second 384-well plate.

**Table 13** Position of ECD digestion reactions for obtaining validation samples

Restriction Enzyme Master Mix to be Validated	Position of ECD Digestion Reaction in 384-Well Plates	
	1-6 Column Runs	12-Column Runs (two 384-well plates produced)
RE Master Mix A	A1	A1 (plate 1)
RE Master Mix B	A2	A2 (plate 1)
RE Master Mix C	B1	B1 (plate 1)
RE Master Mix D	B2	B2 (plate 1)
RE Master Mix E	A13	A1 (plate 2)
RE Master Mix F	A14	A2 (plate 2)
RE Master Mix G	B13	B1 (plate 2)
RE Master Mix H	B14	B2 (plate 2)

- b** Incubate the removed 4- $\mu$ L samples at 80°C for 5 minutes to inactivate the restriction enzymes.
- c** Add 4  $\mu$ L nuclease-free water to each ECD digestion reaction well to restore 10  $\mu$ L/well reaction volumes for the subsequent workflow steps.
- d** Analyze the prepared samples using microfluidic electrophoresis on the 2100 Bioanalyzer (see [page 40](#)) or on the 2200 TapeStation (see [page 41](#)) or using gel electrophoresis (see [page 42](#)).

#### NOTE

After validation of the restriction digest protocol, the ECD sample (from well A1 of the starting 96-well DNA sample plate) is not used for subsequent steps in the workflow.

## Step 1. Digest genomic DNA with restriction enzymes

The ECD sample contains genomic DNA mixed with an 800-bp PCR product that contains restriction sites for all the enzymes used in the digestion protocol. When analyzing validation results, the undigested control should have gDNA bands at >2.5 kbp and a PCR product band at 800 bp. Each of the eight digested ECD samples should have a smear of gDNA restriction fragments between 100 and 2500 bp, overlaid with three predominant bands at approximately 125, 225 and 450 bp. These three bands correspond to the 800-bp PCR product-derived restriction fragments, and precise sizes will differ after digestion in each of the eight RE master mixes.

**NOTE**

In addition to the three predominant bands at approximately 125, 225 and 450 bp, you may detect additional, minor bands in the digested ECD sample lanes.

Successful digestion is indicated by the appearance of the three predominant bands. The presence of additional minor bands, with relative abundance similar to the additional bands visible in [Figure 4](#), [Figure 5](#) and [Figure 6](#), does not impact enrichment results.

It is acceptable for band intensities in digestion reaction B to be slightly reduced, compared to the other digestion reactions.

---

### 3 Sample Preparation

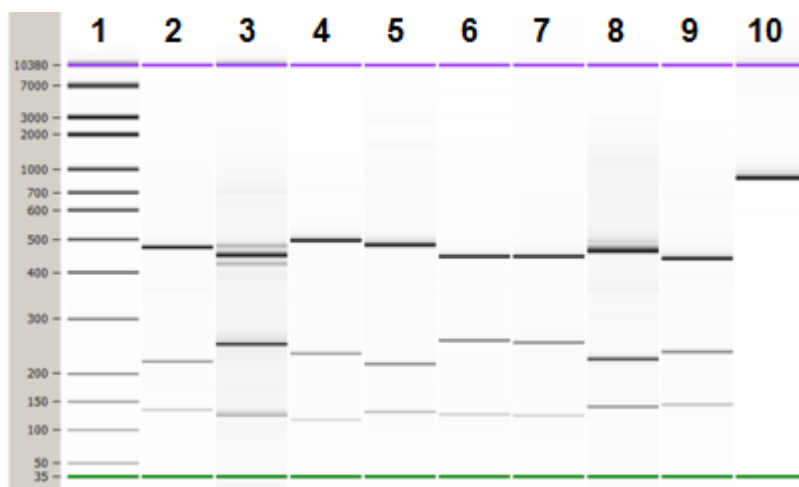
#### Step 1. Digest genomic DNA with restriction enzymes

##### Option 1: Validation by 2100 Bioanalyzer analysis

Use a High Sensitivity DNA Kit (p/n 5067-4626) and the 2100 Bioanalyzer system with 2100 Expert Software (version B.02.07 or higher required to run the High Sensitivity Kit). See the reagent kit guide for general Bioanalyzer system setup instructions.

- Prepare an undigested DNA gel control by combining 0.5  $\mu\text{L}$  of the Enrichment Control DNA stock solution and 3.5  $\mu\text{L}$  of nuclease-free water.
- Prepare the chip, samples and ladder as instructed in the reagent kit guide, using 1  $\mu\text{L}$  of each ECD sample and undigested DNA control for the analysis.
- When loading samples on the chip, load the DNA ladder in the ladder sample well marked on the chip. Load the eight ECD digest samples (A to H) in sample wells 1 to 8, and load the undigested ECD sample in sample well 9. Do not run the undigested ECD control in sample well 1.
- Place the prepared chip into the 2100 Bioanalyzer instrument and start the run within five minutes after preparation.

See [Figure 4](#) for sample Bioanalyzer electrophoresis results.



**Figure 4** Validation of restriction digestion by 2100 Bioanalyzer system analysis. Lane 1: 50-bp DNA ladder, Lanes 2-9: ECD digestion reactions A–H, Lane 10: Undigested Enrichment Control DNA.



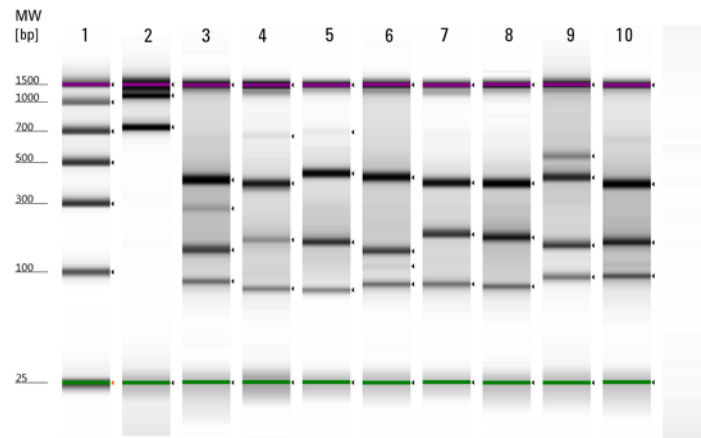
## Step 1. Digest genomic DNA with restriction enzymes

**Option 2: Validation by 2200 TapeStation analysis**

Use a High Sensitivity D1000 ScreenTape (p/n 5067-5584) and reagent kit (p/n 5067-5585). For more information to do this step, see the *2200 TapeStation User Manual*.

- Prepare an undigested DNA gel control by combining 1  $\mu$ L of the Enrichment Control DNA stock solution and 1  $\mu$ L of nuclease-free water.
- Prepare the TapeStation samples as instructed in the *2200 TapeStation User Manual*. Use 2  $\mu$ L of each ECD sample diluted with 2  $\mu$ L of High Sensitivity D1000 sample buffer in separate wells of a tube strip for the analysis.
- Load the sample tube strip, the High Sensitivity D1000 ScreenTape, and loading tips into the 2200 TapeStation as instructed in the *2200 TapeStation User Manual*. Start the run.

See [Figure 5](#) for sample TapeStation electrophoresis results.



**Figure 5** Validation of restriction digestion by 2200 TapeStation analysis. Lane 1: TapeStation D1000 High-Sensitivity Ladder, Lane 2: Undigested Enrichment Control DNA, Lanes 3–10: ECD digestion reactions A–H.

### 3 Sample Preparation

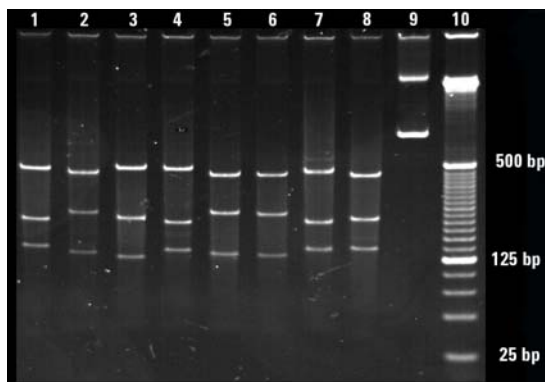
#### Step 1. Digest genomic DNA with restriction enzymes

##### Option 3: Validation by gel electrophoresis

Use a Novex 6% polyacrylamide TBE pre-cast gel and 1X Novex TBE Running Buffer. For more information to do this step, consult the manufacturer's recommendations.

- Prepare an undigested DNA gel control by combining 2  $\mu$ L of the Enrichment Control DNA stock solution and 2  $\mu$ L of nuclease-free water.
- Add 1  $\mu$ L of Novex Hi-Density TBE Sample Buffer (5X) to each 4- $\mu$ L ECD sample.
- Load 5  $\mu$ L of each sample on the gel. In one or more adjacent lanes, load 200 ng of a 50-bp DNA ladder.
- Run the gel at 210 V for approximately 15 minutes.
- Stain the gel in 3X GelRed Nucleic Acid Stain for 10 minutes, and visualize bands under UV radiation.

See [Figure 6](#) for sample gel results.



**Figure 6** Validation of restriction digestion by gel electrophoresis. Lanes 1–8: ECD digestion reactions A–H, Lane 9: Undigested Enrichment Control DNA, Lane 10: 25-bp DNA ladder.

**Stopping Point** If you do not continue to the next step, samples may be stored at  $-20^{\circ}\text{C}$  for long term storage. There are no more long-term stopping points until after the PCR amplification step on [page 72](#).

## Step 2. Hybridize HaloPlex probes A-H to DNA digests A-H for target enrichment and sample indexing

The HaloPlex Exome probe is provided as eight separate probe solutions in wells A-H of the HaloPlex Probe 8-well Strip.

### CAUTION

Do **not** pool the probe solutions provided in the HaloPlex Probe 8-well Strip. The eight probe pools must be used for hybridization in separate wells, with the captured DNA being combined after hybridization.

---

In this step, the eight probe solutions A through H are hybridized, in eight separate hybridization wells, to the DNA restriction fragment libraries A-H prepared in Step 1. The automation protocol below does the liquid handling steps to prepare hybridization reactions in which digests A-H are matched with corresponding probes A-H. (Sample digest A is hybridized to Probe A, while sample digest B is hybridized to Probe B, and so on.)

Hybridization reactions are set up to include the appropriate indexing primers. During the hybridization reaction Illumina sequencing motifs, including index sequences, are integrated into the targeted fragments. For sample indexing primer assignments, see the [Reference](#) chapter for the nucleotide sequences of the 96 indexes used in the HaloPlex Exome Target Enrichment System.

### CAUTION

This guide includes information for kits containing two different sets of indexing primers. **Verify that you are referencing the information appropriate for your kit version before you proceed.**

Kits with indexing primers supplied in a blue plate include 8-bp indexes A01 through H12. See [page 94](#) through [page 95](#) for indexing primer A01–H12 plate map and nucleotide sequence information.

Kits with indexing primers supplied in a clear plate include 8-bp indexes 1 through 96. See [page 98](#) through [page 104](#) for indexing primer 1–96 plate map and nucleotide sequence information.

Protocol steps for indexing using primers provided in either configuration are identical.

---

### 3 Sample Preparation

#### Step 2. Hybridize HaloPlex probes A-H to DNA digests A-H for target enrichment and sample indexing

For 12-column runs, restriction digests from each of the two 384-well plates prepared in Step 1 will be processed during separate segments of the Hybridization\_Exome.pro protocol. Begin the protocol with the first 384-well restriction digest plate on the Bravo deck. The VWorks software will prompt you to replace the first 384-well restriction digest plate with the second restriction digest plate at the appropriate time.

#### Prepare the NGS Bravo

- 1 Gently wipe down the Bravo deck with a NucleoClean decontamination wipe.
- 2 Place a red insert on Bravo deck position 9.
- 3 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.
- 4 Place a 384-well adapter insert on Bravo deck positions 4 and 6.
- 5 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](#).

#### NOTE

To expedite thermal cycler warm-up for the hybridization reaction on [page 50](#), you can enter and initiate the hybridization program on the thermal cycler now, and then pause the program until you are instructed to transfer the reaction plate. Be sure to pause the thermal cycler before the initiation of the 95°C incubation segment. Release the pause immediately after transferring the plate to the thermal cycler for hybridization. Be sure that the 96-well block is in the thermal cycler before initiating the program for warm-up.

## Step 2. Hybridize HaloPlex probes A-H to DNA digests A-H for target enrichment and sample indexing

**Prepare the Indexing Primer Cassette Source Plate****CAUTION**

This guide includes information for kits containing two different sets of indexing primers. **Verify that you are referencing the information appropriate for your kit version before you proceed.**

Reference information for kits with indexing primers supplied in a blue plate begins on [page 92](#).

Reference information for kits with indexing primers supplied in a clear plate begins on [page 96](#).

Protocol steps for indexing using primers provided in either configuration are identical.

---

The Indexing Primer Cassette solutions supplied with the kit must be diluted for use in the exome automation protocol as described below. The well position for each index should correspond to the position of the DNA sample assigned to that index in the original, 96-well pre-digestion DNA sample plate.

- 1 In a full-skirted 96-well Eppendorf twin.tec plate, dispense 22.5  $\mu$ L of nuclease-free water in each sample well used in the run.
- 2 Dispense 27.5  $\mu$ L of the appropriate Indexing Primer Cassette to each sample well position.

Be sure to add only one specific Indexing Primer Cassette to each well, using different indexes for each sample to be multiplexed. Record the identity of the Indexing Primer Cassette assigned to each well for later sequence analysis.

- 3 Seal the plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 3.0 sec. Vortex the plate to mix the index dilutions, then briefly spin to collect the liquid. Keep the Indexing Primer Cassette source plate on ice.

### 3 Sample Preparation

#### Step 2. Hybridize HaloPlex probes A-H to DNA digests A-H for target enrichment and sample indexing

##### Prepare the Hybridization Master Mix Source Plate

- 1 Obtain the black-marked Exome Probe Strip from Box 1. Label the black-marked tube with A, then continue labeling the remaining tubes with B through H, in order. Keep the probe strip tube on ice.

#### CAUTION

It is important to use the probe strip in the proper orientation when preparing the Hybridization Master Mixes as described below. The black color marker on the tube strip and cap strip are used to mark tube A.

- 2 To prepare the eight Hybridization Master Mixes, label eight 1.5-mL tubes with A through H. Dispense the appropriate volume of Hybridization Solution for your run size, shown in [Table 14](#), to each of the eight tubes.

**Table 14** Hybridization Solution volumes for Hybridization Master Mix preparation

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
Hybridization Solution	6.25 µL	79.7 µL	132.8 µL	185.9 µL	239.1 µL	345.3 µL	690.6 µL

- 3 Add the appropriate volume of probe solution from the HaloPlex Probe Strip to the corresponding Hybridization Master Mix tube. Add probe from tube A of the probe strip only to Hybridization Master Mix tube A. Repeat the process for probes B through H, adding each probe solution to a single master mix tube.

**Table 15** HaloPlex Probe volumes for Hybridization Master Mix preparation

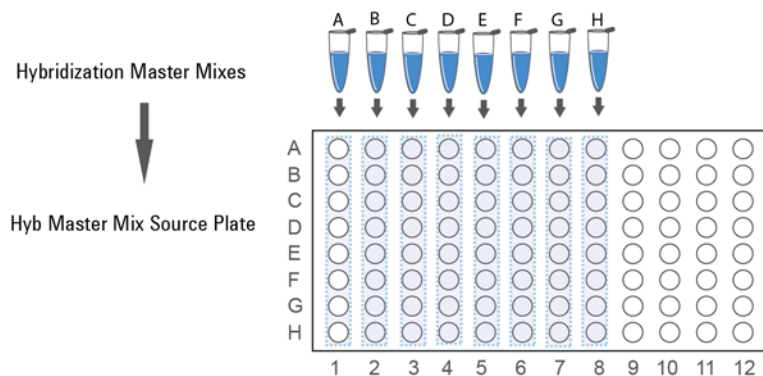
Reagent	Volume for 1 Library	Volume for 1 Column	Volume for 2 Columns	Volume for 3 Columns	Volume for 4 Columns	Volume for 6 Columns	Volume for 12 Columns
HaloPlex Probe A, B, C, D, E, F, G, <b>OR</b> H from provided strip tube	1 µL	12.8 µL	21.3 µL	29.8 µL	38.3 µL	55.3 µL	110.5 µL

## Step 2. Hybridize HaloPlex probes A-H to DNA digests A-H for target enrichment and sample indexing

- 4 Aliquot the Hybridization Master Mixes A through H to a full-skirted 96-well Eppendorf twin.tec plate as shown in Figure 7. Add the volumes indicated in Table 16 of each master mix to each well of the indicated column of the twin.tec plate. Keep the master mixes on ice during the aliquoting steps.

**Table 16** Preparation of the Hybridization Master Mix Source Plate for Hybridization\_Exome.pro

Master Mix Solution	Position on Source Plate	Volume of Master Mix added per Well of Source Plate					
		1-Column Runs	2-Column Runs	3-Column Runs	4-Column Runs	6-Column Runs	12-Column Runs
Hyb Master Mix A	Column 1 (A1-H1)	10.6 $\mu$ L	18.4 $\mu$ L	26.1 $\mu$ L	33.8 $\mu$ L	49.2 $\mu$ L	99.2 $\mu$ L
Hyb Master Mix B	Column 2 (A2-H2)	10.6 $\mu$ L	18.4 $\mu$ L	26.1 $\mu$ L	33.8 $\mu$ L	49.2 $\mu$ L	99.2 $\mu$ L
Hyb Master Mix C	Column 3 (A3-H3)	10.6 $\mu$ L	18.4 $\mu$ L	26.1 $\mu$ L	33.8 $\mu$ L	49.2 $\mu$ L	99.2 $\mu$ L
Hyb Master Mix D	Column 4 (A4-H4)	10.6 $\mu$ L	18.4 $\mu$ L	26.1 $\mu$ L	33.8 $\mu$ L	49.2 $\mu$ L	99.2 $\mu$ L
Hyb Master Mix E	Column 5 (A5-H5)	10.6 $\mu$ L	18.4 $\mu$ L	26.1 $\mu$ L	33.8 $\mu$ L	49.2 $\mu$ L	99.2 $\mu$ L
Hyb Master Mix F	Column 6 (A6-H6)	10.6 $\mu$ L	18.4 $\mu$ L	26.1 $\mu$ L	33.8 $\mu$ L	49.2 $\mu$ L	99.2 $\mu$ L
Hyb Master Mix G	Column 7 (A7-H7)	10.6 $\mu$ L	18.4 $\mu$ L	26.1 $\mu$ L	33.8 $\mu$ L	49.2 $\mu$ L	99.2 $\mu$ L
Hyb Master Mix H	Column 8 (A8-H8)	10.6 $\mu$ L	18.4 $\mu$ L	26.1 $\mu$ L	33.8 $\mu$ L	49.2 $\mu$ L	99.2 $\mu$ L



**Figure 7** Preparation of the Hybridization Master Mix source plate for automation protocol Hybridization\_Exome.pro.

### 3 Sample Preparation

#### Step 2. Hybridize HaloPlex probes A-H to DNA digests A-H for target enrichment and sample indexing

##### Load the NGS Bravo and Run the Hybridization\_Exome.pro VWorks Protocol

- 1 On the VWorks HaloPlex form, under **Step**, select **02 Hybridization\_Exome.pro**.
- 2 Select the number of columns of samples to be processed. Runs must include 1, 2, 3, 4, 6, or 12 columns.
- 3 Click **Update layout and information**.
- 4 Load the Bravo deck according to [Table 17](#).

**Table 17** Initial Bravo deck configuration for Hybridization\_Exome.pro

Location	Content
1	—(empty)—
2	New tip box
3	For 12-column runs only: Empty 384-well Eppendorf twin.tec plate (no 384-well insert required) For 1- to 6-column runs: empty
4	Digested DNA in 384-well plate, seated on 384-well insert *
5	—(empty)—
6	Empty 384-well Eppendorf twin.tec plate, seated on 384-well insert
7	—(empty)—
8	Empty tip box
9	Hybridization Master Mix source plate (96-well Eppendorf twin.tec plate) seated on red insert

\* For 12-column runs, begin the protocol with the first digested DNA sample plate at position 4, and keep the second digested DNA sample plate sealed on ice until you are prompted to add it to the NGS Bravo during the run.

- 5 Verify that the NGS Bravo has been set up as displayed in the **Bravo Deck Setup** and **Information** regions of the form.
- 6 Verify that the **Current Tip State** indicator on the form matches the configuration of unused and used tips in the tip boxes at Bravo Deck positions 2 and 8, respectively. See [page 19](#) for more information on using this segment of the form during the run.
- 7 When verification is complete, click **Start** to start the run.



## Step 2. Hybridize HaloPlex probes A-H to DNA digests A-H for target enrichment and sample indexing

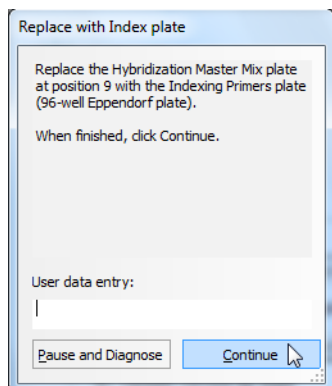


For each DNA sample, the NGS Bravo combines the completed RE digestion reaction A-H with the corresponding Hybridization Master Mix A-H in wells of a 384-well plate.

The remaining steps of the **Hybridization\_Exome.pro** automation protocol differ for runs of 1-6 columns of samples (where a single 384-well digested DNA sample plate is processed) versus runs of 12 columns (where two 384-well digested DNA sample plates are processed). For 1-6 column runs, follow the steps immediately below. For 12-column runs, go to [page 51](#).

### Hybridization Workflow for 1-6 Column Runs

- 1 When the NGS Bravo has finished combining the DNA digests and Hybridization master mixes, you will be prompted by VWorks as shown below.

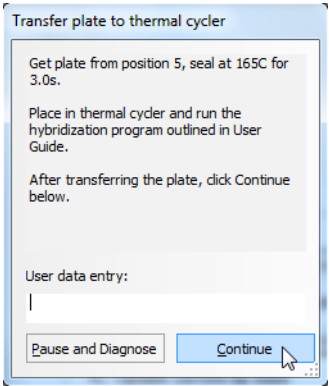


- a Remove and discard the Hybridization Master Mix source plate from position 9.
- b Obtain the Indexing Primer Cassette source plate (prepared on [page 45](#)), carefully unseal the plate to prevent splashing, and load the plate on position 9, seated on the red insert.
- c When finished, click **Continue** on the VWorks prompt.

### 3 Sample Preparation

#### Step 2. Hybridize HaloPlex probes A-H to DNA digests A-H for target enrichment and sample indexing

- 2 When the NGS Bravo has finished adding the indexing primers to the hybridization reactions, you will be prompted by VWorks as shown below.



- a Remove the hybridization plate from position 5. Seal the sample plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 3.0 sec. Spin the plate briefly.
- b When finished, click **Continue** on the VWorks prompt.
- c Place the plate in the thermal cycler and run the hybridization program in [Table 18](#). Use a heated lid. Do **not** include a low-temperature hold step in the thermal cycler program. Incubation at 54°C for greater than 24 hours is not recommended.

**Table 18** Thermal cycler program for HaloPlex Exome probe hybridization

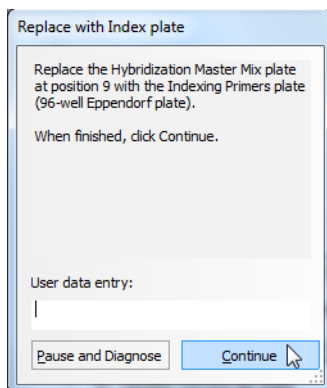
Step	Temperature	Time
Step 1	95°C	10 minutes
Step 2	54°C	16-24 hours

- d After the 16-24 hour incubation step, proceed to “[Step 3. Pool hybridized DNA samples](#)” on page 55.

## Step 2. Hybridize HaloPlex probes A-H to DNA digests A-H for target enrichment and sample indexing

**Hybridization Workflow for 12 Column Runs**

- 1 When the NGS Bravo has finished combining the DNA digests and Hybridization master mixes for the first plate of digested DNA samples, you will be prompted by VWorks as shown below.

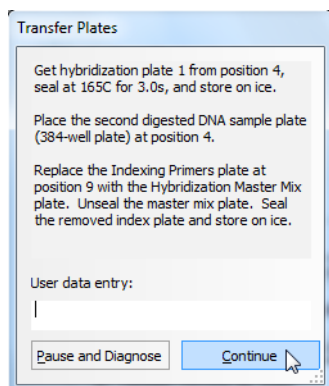


- a Remove the Hybridization Master Mix source plate from position 9. Seal the master mix plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 3.0 sec, and store the plate on ice.
- b Obtain the Indexing Primer source plate (prepared on [page 45](#)), carefully unseal the plate to prevent splashing, and load the plate on position 9, seated on the red insert.
- c When finished, click **Continue** on the VWorks prompt.

### 3 Sample Preparation

#### Step 2. Hybridize HaloPlex probes A-H to DNA digests A-H for target enrichment and sample indexing

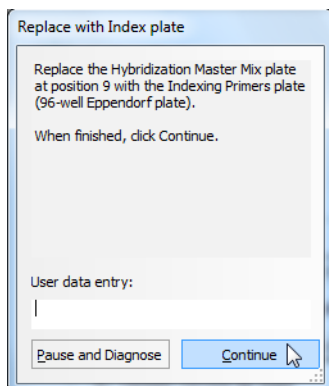
- 2 When the NGS Bravo has finished adding the indexing primers to the first plate of hybridization reactions, you will be prompted by VWorks as shown below.



- a Remove the hybridization plate 1 from position 4. Seal the plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 3.0 sec. Place hybridization plate 1 on ice while the NGS Bravo prepares the second hybridization plate.
- b Place the second digested DNA sample plate at position 4, seated on the 384-well plate insert. Carefully unseal the plate.
- c Remove the Indexing Primer source plate from position 9. Seal the plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 3.0 sec, and store the plate on ice.
- d Obtain the Hybridization Master Mix source plate from ice, carefully unseal the plate, and load the plate on position 9, seated on the red insert.
- e When finished, click **Continue** on the VWorks prompt.

## Step 2. Hybridize HaloPlex probes A-H to DNA digests A-H for target enrichment and sample indexing

- 3 When the NGS Bravo has finished combining the DNA digests and Hybridization master mixes for the second plate of digested DNA samples, you will be prompted by VWorks as shown below.

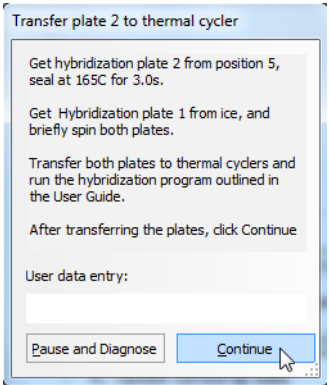


- a Remove and discard the Hybridization Master Mix source plate from position 9.
- b Obtain the Indexing Primer source plate from storage on ice, carefully unseal the plate, and load the plate on position 9, seated on the red insert.
- c When finished, click **Continue** on the VWorks prompt.

### 3 Sample Preparation

#### Step 2. Hybridize HaloPlex probes A-H to DNA digests A-H for target enrichment and sample indexing

- 4 When the NGS Bravo has finished adding the indexing primers to the second plate of hybridization reactions, you will be prompted by VWorks as shown below.



- a Remove the second hybridization plate from position 5. Seal the plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 3.0 sec.
- b Recover the first hybridization plate from ice storage. Spin both hybridization plates briefly, then place each plate in a thermal cycler and run the hybridization program in [Table 19](#). Use a heated lid. Do **not** include a low-temperature hold step in the thermal cycler program. Incubation at 54°C for greater than 24 hours is not recommended.

**Table 19** Thermal cycler program for HaloPlex Exome probe hybridization

Step	Temperature	Time
Step 1	95°C	10 minutes
Step 2	54°C	16–24 hours

- c When finished transferring the plates, click **Continue** on the VWorks prompt.
- d After completing the 16-24 hour incubation, proceed to “[Step 3. Pool hybridized DNA samples](#)” on page 55.

## Step 3. Pool hybridized DNA samples

In this step, the eight hybridization reactions that correspond to one DNA sample are combined into a single well and mixed with streptavidin beads in preparation for the capture protocol.

The NGS Bravo pools the hybridized DNA samples from the 384-well hybridization plates into wells of a Nunc DeepWell plate in the same sample configuration that was used for the pre-digestion DNA sample plate.

### Prepare the NGS Bravo and reagents

- 1** Remove reagents to be used in this protocol from cold storage and allow the solutions to reach room temperature:
  - From  $-20^{\circ}\text{C}$  storage, remove the Capture Solution.
  - From  $+4^{\circ}\text{C}$  storage, remove the HaloPlex Magnetic Beads.
- 2** Gently wipe down the Bravo deck with a NucleoClean decontamination wipe.
- 3** For all run sizes, place a 384-well adapter insert on Bravo deck position 6. For 12-column runs only, place a second 384-well adapter insert on Bravo deck position 4.
- 4** Pre-set the temperature of Bravo deck positions 4 and 6 to  $54^{\circ}\text{C}$  using the Inheco Multi TEC control touchscreen, as described in [Setting the Temperature of Bravo Deck Heat Blocks](#).

**3 Sample Preparation**  
Step 3. Pool hybridized DNA samples

**Prepare the HaloPlex Magnetic Beads Source Plate**

- 1** Vigorously resuspend the HaloPlex Magnetic Beads on a vortex mixer. The beads settle during storage.
- 2** Wash the magnetic beads.
  - a** Transfer 40  $\mu$ L per sample of the HaloPlex Magnetic Beads suspension to a 1.5-mL tube or conical vial, using volumes provided in [Table 20](#).

**Table 20** Volume of HaloPlex Magnetic Bead suspension for Post Hybridization Pooling\_Exome.pro

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
HaloPlex Magnetic Beads	0.04 mL	0.36 mL	0.68 mL	1.0 mL	1.32 mL	1.96 mL	3.92 mL

- b** Put the vial into a compatible magnetic device for 5 minutes.
- c** After verifying that the solution has cleared, carefully remove and discard the supernatant using a pipette.
- d** Add an equivalent volume of Capture Solution (see [Table 21](#)) to the beads and resuspend by pipetting up and down.

**Table 21** Volume of Capture Solution used for bead resuspension

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
Capture Solution	0.04 mL	0.36 mL	0.68 mL	1.0 mL	1.32 mL	1.96 mL	3.92 mL

- 3** Prepare a Nunc DeepWell source plate for the washed HaloPlex streptavidin bead suspension. Add 40  $\mu$ L of the homogeneous bead suspension to all wells of the Nunc DeepWell plate that correspond to sample-containing wells of the pre-digestion DNA sample plate.



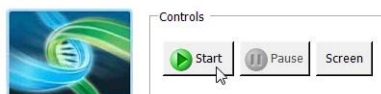
### Load the NGS Bravo and Run the Post Hybridization Pooling\_Exome.pro VWorks Protocol

- 1 On the VWorks HaloPlex form, under **Step**, select **03 Post Hybridization Pooling\_Exome.pro**.
- 2 Select the number of columns of samples to be processed. Runs must include 1, 2, 3, 4, 6, or 12 columns.
- 3 Click **Update layout and information**.
- 4 Load the Bravo deck according to [Table 22](#).

**Table 22** Initial Bravo deck configuration for Post Hybridization Pooling\_Exome.pro

Location	Content
1	—(empty)—
2	New tip box
3	—(empty)—
4	384-well adapter insert
5	HaloPlex magnetic streptavidin bead plate (Nunc DeepWell plate)
6	384-well adapter insert
7	—(empty)—
8	Empty tip box
9	—(empty)—

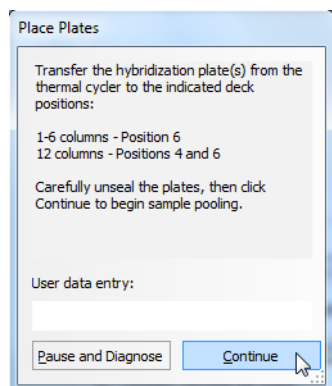
- 5 Verify that the NGS Bravo has been set up as displayed in the **Bravo Deck Setup** and **Information** regions of the form.
- 6 Verify that the **Current Tip State** indicator on the form matches the configuration of unused and used tips in the tip boxes at Bravo Deck positions 2 and 8, respectively. See [page 19](#) for more information on using this segment of the form during the run.
- 7 When verification is complete, click **Start** to start the run.



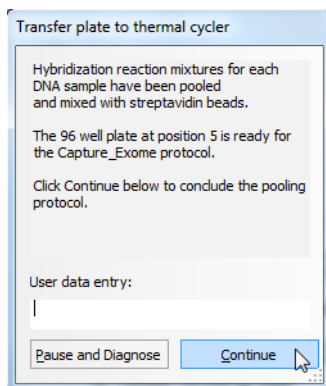
### 3 Sample Preparation

#### Step 3. Pool hybridized DNA samples

- 8 When prompted by VWorks as shown below, get the hybridization plate(s) from the thermal cycler. Briefly spin the sealed plates to collect the liquid, then load the plate(s) on the indicated position(s) of the Bravo deck. Carefully unseal the plate(s). When finished, Click **Continue** on the prompt to resume the pooling protocol.



The NGS Bravo pools the target DNA-HaloPlex probe hybrids into the appropriate wells of the streptavidin bead-containing Nunc DeepWell plate. When the Bravo has finished the pooling step, you will be prompted by VWorks as shown below.



Remove the sample capture plate from position 5, seal the plate, and store at room temperature while you prepare the components of the Capture\_Exome automation protocol. Click **Continue** on the prompt to conclude the pooling protocol.

## Step 4. Capture and amplify the target DNA

In this step, the pooled target DNA-HaloPlex probe hybrids are captured on streptavidin beads. After capture, DNA ligase is added to seal nicks, then target DNA is eluted and PCR-amplified.

### Assemble reagents for the run

- 1 Remove the Wash Solution, Ligation Solution and SSC Buffer from  $-20^{\circ}\text{C}$  storage and allow the solutions to reach room temperature.
- 2 Prepare 50  $\mu\text{L}$  per sample, plus excess, of fresh 50 mM NaOH for use in the DNA elution step on [page 60](#).

Prepare the 50 mM NaOH solution from a 10M NaOH stock solution.

### CAUTION

Using high-quality NaOH is critical for optimal DNA elution and recovery.

- Do not use stock NaOH solutions that were stored at concentrations below 10 M to prepare the 50 mM NaOH solution.
- Keep the 50 mM NaOH solution container sealed when not in use, especially when processing large numbers of samples per run.

**Table 23** Amount of 50mM NaOH required per run size

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
60 $\mu\text{L}$	540 $\mu\text{L}$	1020 $\mu\text{L}$	1500 $\mu\text{L}$	1980 $\mu\text{L}$	2940 $\mu\text{L}$	5880 $\mu\text{L}$

- 3 Obtain or prepare 1  $\mu\text{L}$  per sample, plus excess, of 2 M acetic acid, for use in the PCR master mix on [page 61](#).

### CAUTION

It is critical to use high-quality acetic acid at 2 M concentration in this step to ensure neutralization of the NaOH used for elution.

See [Table 1](#) on page 9 for 2 M acetic acid supplier information, or prepare 2 M acetic acid from high-quality glacial acetic acid.

### 3 Sample Preparation

#### Step 4. Capture and amplify the target DNA

##### Prepare the NGS Bravo

- 1 Gently wipe down the Bravo deck with a NucleoClean decontamination wipe.
- 2 Place a red insert on Bravo deck position 4.
- 3 Pre-set the temperature of Bravo deck position 4 to 46°C using the Inheco Multi TEC control touchscreen, as described in [Setting the Temperature of Bravo Deck Heat Blocks](#).
- 4 Place a second red insert on Bravo deck position 6.
- 5 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](#).
- 6 Place the silver Nunc plate insert on Bravo deck position 9.
- 7 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.

##### Prepare wash and elution solution source plates

Prepare a separate source plate for each of the solutions listed in [Table 24](#). Use full-skirted 96-well Eppendorf twin.tec plates to prepare all three source plates. For all sample-containing well positions, add the specified volume of solution to each source plate.

**Table 24** Preparation of solution source plates for Capture\_Exome.pro protocol

Solution	Volume to dispense per well of source plate
Wash Solution	110 µL
SSC Buffer	110 µL
50 mM NaOH	60 µL

Seal the 50 mM NaOH source plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 1.0 sec. Leave sealed until you are prompted to add the plate to the Bravo deck in [step 14](#) on [page 70](#).

**Prepare the Master Mixes for Capture\_Exome.pro protocol**

- 1 Prepare the appropriate amount of PCR Master Mix, according to the table below.

Mix well by gentle vortexing, then spin the tube briefly.

**Table 25** Preparation of PCR Master Mix for Capture\_Exome.pro

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	32.2 µL	410.6 µL	684.3 µL	958.0 µL	1231.7 µL	1779.1 µL	3558.1 µL
5X Herculase II Reaction Buffer	20 µL	255.0 µL	425.0 µL	595.0 µL	765.0 µL	1105.0 µL	2210.0 µL
dNTPs (100 mM)	0.8 µL	10.2 µL	17.0 µL	23.8 µL	30.6 µL	44.2 µL	88.4 µL
Primer 1	2 µL	25.5 µL	42.5 µL	59.5 µL	76.5 µL	110.5 µL	221.0 µL
Primer 2	2 µL	25.5 µL	42.5 µL	59.5 µL	76.5 µL	110.5 µL	221.0 µL
2 M Acetic acid	1 µL	12.8 µL	21.3 µL	29.8 µL	38.3 µL	55.3 µL	110.5 µL
Herculase II Fusion DNA Polymerase	2 µL	25.5 µL	42.5 µL	59.5 µL	76.5 µL	110.5 µL	221.0 µL
<b>Total Volume</b>	<b>60 µL</b>	<b>765 µL</b>	<b>1275 µL</b>	<b>1785 µL</b>	<b>2295 µL</b>	<b>3315 µL</b>	<b>6630 µL</b>

### 3 Sample Preparation

#### Step 4. Capture and amplify the target DNA

- 2 Prepare the appropriate amount of Ligation Master Mix, according to the table below.

Mix well by gentle vortexing, then spin the tube briefly. Store the master mix on ice until it is used on [page 66](#). The Ligation Master Mix is added to the Master Mix Source Plate just before it is used in the Capture\_Exome.pro protocol. Do not add this master mix to the source plate before starting the run.

**Table 26** Preparation of Ligation Master Mix for Capture\_Exome.pro

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
Ligation Solution	47.5 µL	605.6 µL	1009 µL	1413 µL	1817 µL	2624 µL	5249 µL
DNA Ligase	2.5 µL	31.9 µL	53.1 µL	74.4 µL	95.6 µL	138.1 µL	276.3 µL
<b>Total Volume</b>	<b>50 µL</b>	<b>637.5 µL</b>	<b>1062.1 µL</b>	<b>1487.4 µL</b>	<b>1912.6 µL</b>	<b>2762.1 µL</b>	<b>5525.3 µL</b>

#### Prepare the Master Mix Source Plate for Capture\_Exome.pro

To prepare the Master Mix source plate for Capture\_Exome.pro, add the volume indicated in [Table 27](#) of PCR Master Mix to all wells of column 3 of a Nunc DeepWell plate.

**Table 27** Preparation of the Master Mix Source Plate for Capture\_Exome.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 Master Mix	Column 3 (A3-H3)	88.1 µL	151.9 µL	215.6 µL	279.4 µL	406.9 µL	821.3 µL

#### NOTE

Columns 1 and 2 of the Master Mix source plate remain empty at this step. You will be prompted to add Ligation Master Mix to Column 2 at the appropriate time during the Capture\_Exome.pro protocol. Column 1 will remain empty during the Capture\_Exome.pro protocol.

**Load the NGS Bravo and Run the Capture\_Exome.pro VWorks Protocol**

- 1 On the VWorks HaloPlex form, under **Step**, select **04 Capture\_Exome.pro**.
- 2 Select the number of columns of samples to be processed. Runs must include 1, 2, 3, 4, 6, or 12 columns.
- 3 Click **Update layout and information**.
- 4 Load the Bravo deck according to [Table 28](#).

**Table 28** Initial Bravo deck configuration for Capture\_Exome.pro

Location	Content
1	Empty Axygen 96 Deep Well Plate (square wells) for waste
2	New tip box
3	Wash Solution source plate (full-skirted 96-well Eppendorf twin.tec plate)
4	Empty half-skirted 96-well Eppendorf twin.tec plate seated on red insert
5	Pooled hybridized DNA samples + streptavidin beads in Nunc DeepWell plate
6	Empty half-skirted 96-well Eppendorf twin.tec plate seated on red insert
7	—(empty)—
8	Empty tip box
9	Master Mix source plate (Nunc DeepWell plate) seated on silver insert

- 5 Verify that the NGS Bravo has been set up as displayed in the **Bravo Deck Setup** and **Information** regions of the form.
- 6 Verify that the **Current Tip State** indicator on the form matches the configuration of unused and used tips in the tip boxes at Bravo Deck positions 2 and 8, respectively. See [page 19](#) for more information on using this segment of the form during the run.
- 7 When verification is complete, click **Start** to start the run.



### 3 Sample Preparation

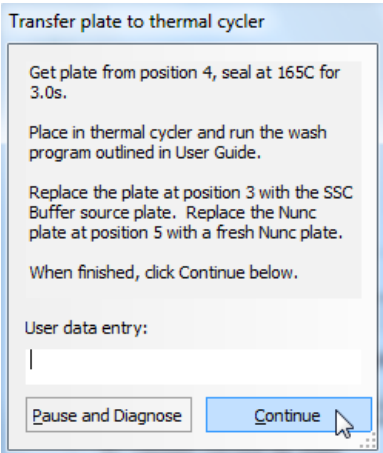
#### Step 4. Capture and amplify the target DNA

The NGS Bravo completes the liquid-handling steps for capture and washing of the target DNA-HaloPlex probe hybrids on the streptavidin beads.

#### NOTE

To expedite thermal cycler warm-up for the subsequent wash program on [page 64](#), you can enter and initiate the wash program on the thermal cycler now, and then pause the program until you are instructed to transfer the reaction plate. Be sure to pause the thermal cycler before the initiation of the 46°C incubation segment. Release the pause immediately after transferring the plate to the thermal cycler in [step 8](#).

- 8 When the NGS Bravo has finished preparing the capture wash plate, you will be prompted by VWorks as shown below.



- a Get the sample plate from position 4 and seal the plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 3.0 sec.

Transfer the sealed plate to a thermal cycler and run the wash program shown in [Table 29](#), using a heated lid.

Do **not** include a low-temperature hold step in the thermal cycler program following the 10-minute incubation.

**Table 29** Thermal cycler program for Capture\_Exome.pro wash step

Step	Temperature	Time
Step 1	46°C	10 minutes



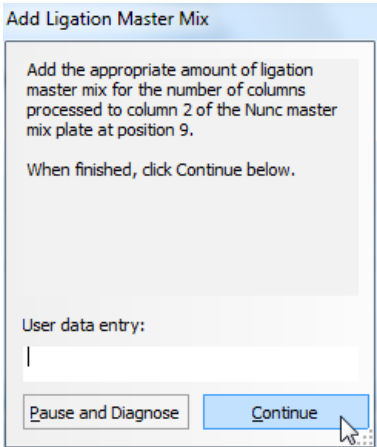
## Step 4. Capture and amplify the target DNA

- b** Remove and discard the Wash Solution plate from position 3 of the Bravo deck. Place the SSC Buffer source plate at position 3 for use in the ligation segment of the protocol.
- c** Remove and discard the used Nunc plate from position 5 of the Bravo deck. Place a fresh Nunc DeepWell plate at position 5.
- d** After completing all steps, click **Continue** on the VWorks prompt to continue the automation protocol. Do not wait for the conclusion of the thermal cycler wash program to continue the protocol.

### 3 Sample Preparation

#### Step 4. Capture and amplify the target DNA

- 9 During the 10-minute incubation of the sample plate on the thermal cycler, you will be prompted to add the Ligation Master Mix to the Master Mix source plate as shown below.



Add the volume of Ligation Master Mix indicated in [Table 30](#) to all wells of column 2 of the Nunc DeepWell Master Mix source plate on Bravo deck position 9.

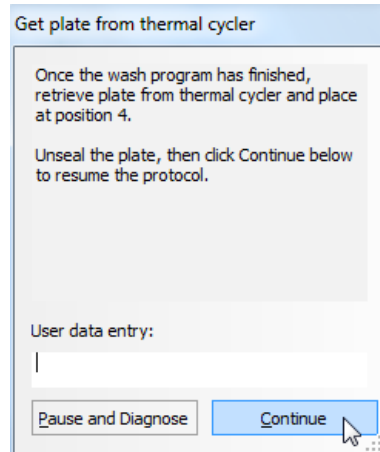
**Table 30** Addition of Ligation Master Mix to the Master Mix Source Plate for Capture\_Exome.pro

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
Ligation Master Mix	Column 2 (A2-H2)	73.4 µL	126.6 µL	179.7 µL	232.8 µL	339.1 µL	684.4 µL

#### NOTE

The Master Mix source plate at position 9 should already contain the PCR Master Mix in Column 3. Be sure to add the Ligation Master Mix to Column 2 of the source plate at this step.

- 10** Once the wash program in [Table 29](#) has finished and you are prompted by VWorks, transfer the plate from the thermal cycler to Bravo deck position 4. Carefully unseal the plate, then click **Continue** on the VWorks prompt to resume the Capture\_Exome.pro protocol.



The NGS Bravo completes the liquid-handling steps for ligation of the captured target DNA.

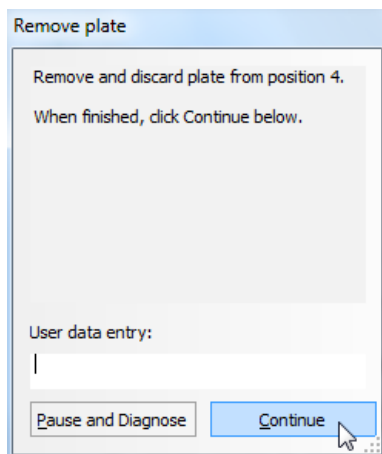
**NOTE**

To expedite thermal cycler warm-up for the subsequent ligation program on [page 69](#), you can enter and initiate the program on the thermal cycler now, and then pause the program until you are instructed to transfer the reaction plate. Be sure to pause the thermal cycler before the initiation of the 55°C incubation segment. Release the pause immediately after transferring the plate to the thermal cycler in [step 12](#).

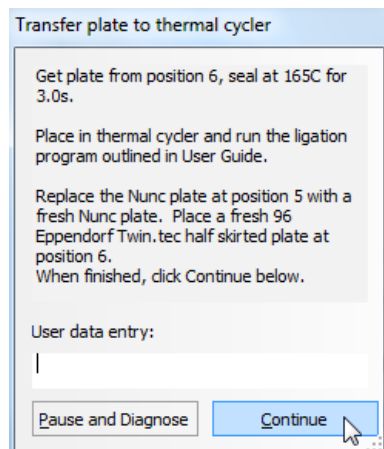
### 3 Sample Preparation

#### Step 4. Capture and amplify the target DNA

- 11 When prompted by VWorks as shown below, remove and discard the plate from position 4. Click **Continue** on the prompt to resume the protocol.



**12** When the NGS Bravo has finished preparing the ligation plate, you will be prompted by VWorks as shown below.



- a** Get the sample plate from position 6 and seal the plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 3.0 sec.
- b** Transfer the sealed plate to a thermal cycler and run the ligation program shown in [Table 31](#), using a heated lid.

**Table 31** Thermal cycler program for Capture\_Exome.pro ligation step

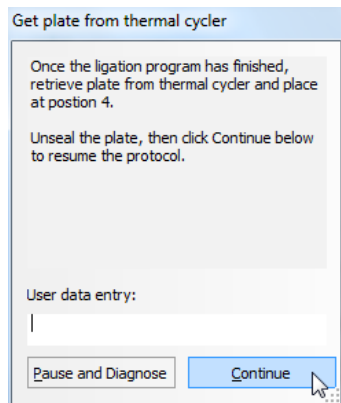
Step	Temperature	Time
Step 1	55°C	10 minutes
Step 2	4°C	Hold

- c** Remove and discard the Nunc DeepWell plate from position 5 of the Bravo deck. Place a fresh Nunc DeepWell plate at position 5 for use in the ligation purification steps of the protocol.
- d** Place a fresh half-skirted 96-well Eppendorf twin.tec plate at position 6 for use in the following PCR protocol steps.
- e** After completing all steps, click **Continue** on the VWorks prompt to continue the automation protocol. Do not wait for the conclusion of the thermal cycler ligation program to continue the protocol.

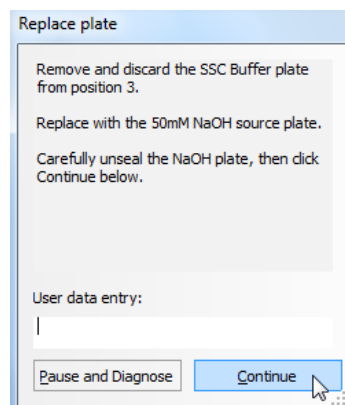
### 3 Sample Preparation

#### Step 4. Capture and amplify the target DNA

- 13** Once the ligation program in [Table 31](#) is finished and you are prompted by VWorks, transfer the plate from the thermal cycler to Bravo deck position 4. Carefully unseal the plate, then click **Continue** on the VWorks prompt to resume the Capture\_Exome.pro protocol.



- 14** When prompted by VWorks as shown below, remove and discard the SSC Buffer plate from position 3 of the Bravo deck. Place the 50 mM NaOH source plate at position 3 for use in the elution steps of the protocol. After carefully unsealing the source plate, click **Continue** on the prompt to resume the protocol.

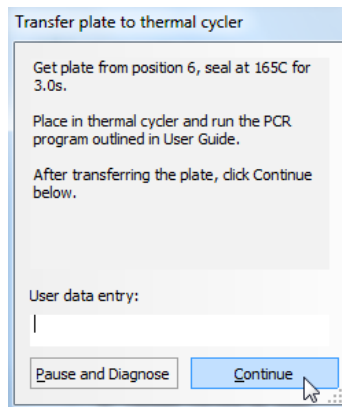


The NGS Bravo completes the liquid-handling steps for elution of the captured target DNA followed by preparation of PCR reactions for amplification.

**NOTE**

To expedite thermal cycler warm-up for the subsequent PCR program on [page 72](#), you can enter and initiate the PCR program on the thermal cycler now, and then pause the program until you are instructed to transfer the reaction plate. Be sure to pause the thermal cycler before the initiation of the 98°C denaturation segment. Release the pause immediately after transferring the plate to the thermal cycler in [step 15](#).

- 15** When the NGS Bravo has finished preparing the PCR amplification reactions, you will be prompted by VWorks as shown below.



- a** Get the sample plate from position 6 and seal the plate using the PlateLoc Thermal Microplate Sealer, with sealing settings of 165°C and 3.0 sec.
- b** Transfer the sealed plate to a thermal cycler and run the PCR program in [Table 32](#), using a heated lid.

### 3 Sample Preparation

#### Step 4. Capture and amplify the target DNA

Consult the Certificate of Analysis (provided with HaloPlex Exome Target Enrichment System Box 1) for the PCR cycling recommendation.

**Table 32** HaloPlex post-capture DNA amplification PCR program

Segment	Number of Cycles	Temperature	Time
1	1	98°C	2 minutes
2	Obtain cycle number from Certificate of Analysis	98°C	30 seconds
		60°C	30 seconds
		72°C	1 minute
3	1	72°C	10 minutes
4	1	8°C	Hold

- c** After initiating the PCR program in the thermal cycler, click **Continue** on the VWorks prompt to finish the automation protocol.
- d** If you are continuing to the next step of PCR product purification, remove the Agencourt AMPure XP Beads from +4°C storage for use on [page 73](#). Let the beads come to room temperature for the remainder of the amplification program.

**Stopping Point** If you do not continue to the next step, PCR products may be stored at –20°C for up to 72 hours or at 8°C overnight. For best results, however, purify PCR products as soon as possible.



## Step 5. Purify the amplified target DNA

In this step, the NGS Bravo does the liquid handling steps to purify the amplified target DNA sample using AMPure XP beads.

### Prepare the NGS Bravo and reagents

- 1 Gently wipe down the Bravo deck with a NucleoClean decontamination wipe.
- 2 Let the AMPure XP beads come to room temperature for at least 30 minutes.  
*Do not freeze the AMPure XP beads at any time.*
- 3 Mix the room-temperature bead suspension well so that the reagent appears homogeneous and consistent in color.
- 4 Prepare a Nunc DeepWell source plate containing AMPure XP beads. For each well to be processed, add 100  $\mu$ L of homogeneous AMPure XP beads per well to the Nunc DeepWell plate.

#### NOTE

After preparing the source plate, retain the beads at room temperature for use in the next step of the workflow on [page 77](#).

- 5 Place a red insert on Bravo deck position 6.
- 6 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](#).
- 7 Prepare a Thermo Scientific reservoir containing 15 mL of nuclease-free water.
- 8 Prepare a Thermo Scientific reservoir containing 15 mL of the final sample elution buffer [nuclease-free 10 mM Tris-acetate or Tris-HCl buffer (pH 8.0)].
- 9 Prepare a separate Thermo Scientific reservoir containing 45 mL of freshly-prepared 70% ethanol.

### Load the NGS Bravo and Run the Purification\_01\_Exome.pro VWorks Protocol

- 1 On the VWorks HaloPlex form, under **Step**, select **05 Purification\_01\_Exome.pro**.
- 2 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 5. Purify the amplified target DNA

- 3 Click **Update layout and information**.
- 4 Load the Bravo deck according to [Table 33](#).

**Table 33** Initial Bravo deck configuration for Purification\_01\_Exome.pro

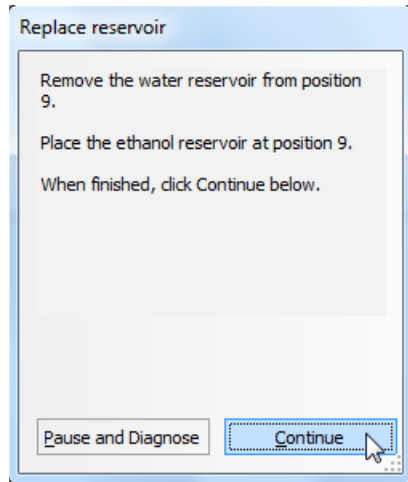
Location	Content
1	Empty Axygen 96 Deep Well Plate (square wells) for waste
2	New tip box
3	—(empty)—
4	—(empty)—
5	AMPure XP beads in Nunc DeepWell source plate
6	Amplified DNA samples in half-skirted 96-well Eppendorf twin.tec plate seated on red insert
7	—(empty)—
8	Empty tip box
9	Nuclease-free water in Thermo Scientific reservoir

- 5 Verify that the NGS Bravo has been set up as displayed in the **Bravo Deck Setup** region of the form.
- 6 Verify that the **Current Tip State** indicator on the form matches the configuration of unused and used tips in the tip boxes at Bravo Deck positions 2 and 8, respectively. See [page 19](#) for more information on using this segment of the form during the run.
- 7 When verification is complete, click **Start** to start the run.



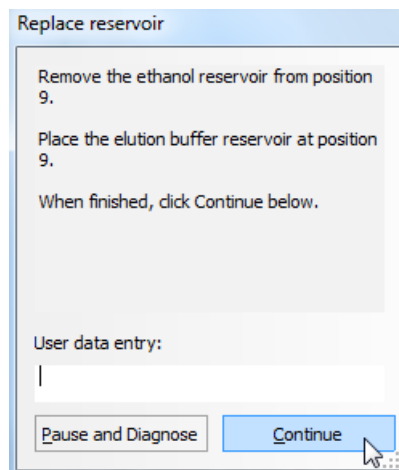
- 8 When prompted by VWorks as shown below, remove the water reservoir from position 9 of the Bravo deck and replace it with the 70% ethanol reservoir.

When finished, click **Continue** on the VWorks prompt.



- 9 When prompted by VWorks as shown below, remove the 70% ethanol reservoir from position 9 of the Bravo deck and replace it with the final sample elution buffer reservoir.

When finished, click **Continue** on the VWorks prompt.



### 3 Sample Preparation

#### Step 5. Purify the amplified target DNA

The NGS Bravo completes the liquid-handling steps for elution of the captured target DNA.

When the automation protocol is finished, the DNA samples are in the bead-containing Nunc DeepWell plate at position 7. Leave the sample plate on the Bravo deck for the subsequent automation protocol,

**Purification\_02\_Exome.pro.**

## Step 6. Remove primer dimers in second serial purification

In this step, the NGS Bravo does the liquid handling steps to remove primer dimers from the enriched DNA library using AMPure XP beads.

- 1 Verify that the AMPure XP beads are at room temperature. If beads are in cold storage, let beads come to room temperature for at least 30 minutes.
- 2 Mix the room-temperature bead suspension well so that the reagent appears homogeneous and consistent in color.
- 3 Prepare a Nunc DeepWell source plate containing AMPure XP beads. For each well to be processed, add 100  $\mu$ L of homogeneous AMPure XP beads per well to the Nunc DeepWell plate.
- 4 Prepare a Thermo Scientific reservoir containing 15 mL of nuclease-free water.
- 5 Prepare a Thermo Scientific reservoir containing 15 mL of the final sample elution buffer [nuclease-free 10 mM Tris-acetate or Tris-HCl buffer (pH 8.0)].
- 6 Prepare a separate Thermo Scientific reservoir containing 45 mL of freshly-prepared 70% ethanol.

### Load the NGS Bravo and Run the Purification\_02\_Exome.pro. VWorks Protocol

- 1 On the VWorks HaloPlex form, under **Step**, select **06 Purification\_02\_Exome.pro**.
- 2 Select the number of columns of samples to be processed. Runs must include 1, 2, 3, 4, 6, or 12 columns.
- 3 Click **Update layout and information**.

### 3 Sample Preparation

#### Step 6. Remove primer dimers in second serial purification

4 Load the Bravo deck according to [Table 34](#).

**Table 34** Initial Bravo deck configuration for Purification\_02\_Exome.pro.

Location	Content
1	Empty Axygen 96 Deep Well Plate (square wells) for waste
2	New tip box
3	Empty full-skirted 96-well Eppendorf twin.tec plate
4	—(empty)—
5	AMPure XP beads in Nunc DeepWell source plate
6	—(empty)—
7	Library samples for purification in Nunc DeepWell plate
8	Empty tip box
9	Nuclease-free water in Thermo Scientific reservoir

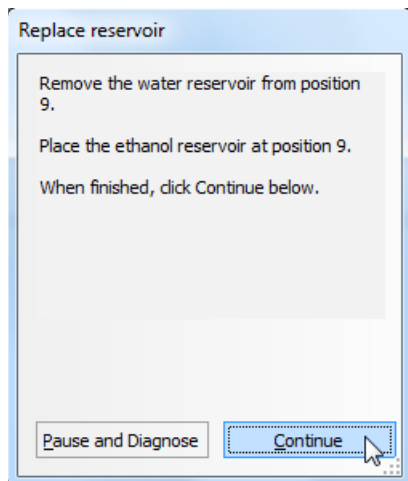
- 5 Verify that the NGS Bravo has been set up as displayed in the **Bravo Deck Setup** region of the form.
- 6 Verify that the **Current Tip State** indicator on the form matches the configuration of unused and used tips in the tip boxes at Bravo Deck positions 2 and 8, respectively. See [page 19](#) for more information on using this segment of the form during the run.
- 7 When verification is complete, click **Start** to start the run.



## Step 6. Remove primer dimers in second serial purification

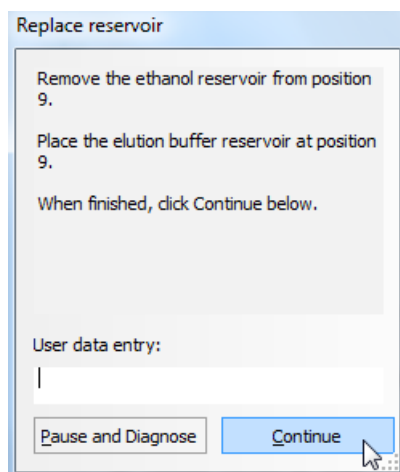
- 8 When prompted by VWorks as shown below, remove the water reservoir from position 9 of the Bravo deck and replace it with the 70% ethanol reservoir.

When finished, click **Continue** on the VWorks prompt.



- 9 When prompted by VWorks as shown below, remove the 70% ethanol reservoir from position 9 of the Bravo deck and replace it with the final sample elution buffer reservoir.

When finished, click **Continue** on the VWorks prompt.

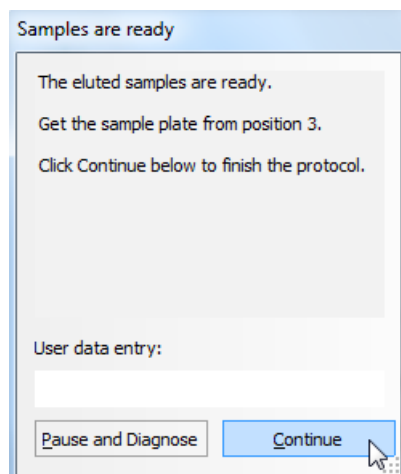


### 3 Sample Preparation

#### Step 6. Remove primer dimers in second serial purification

The NGS Bravo completes the liquid-handling steps for elution of the captured target DNA.

- 10** When the NGS Bravo has finished preparing the final eluted sample plate, you will be prompted by VWorks as shown below. Click **Continue** on the VWorks prompt to finish the protocol.



**Stopping Point** If you do not continue to the next step, samples may be stored at  $-20^{\circ}\text{C}$  for long-term storage (up to one year). Avoid subjecting the stored DNA samples to multiple freeze-thaw cycles.



## Step 7. Validate enrichment and quantify enriched target DNA

Prior to sample pooling and sequencing sample preparation, validate enrichment and quantify the enriched target DNA in each library sample by microfluidics analysis using the 2100 Bioanalyzer (see [page 82](#)) or the 2200 TapeStation (see [page 83](#)).

Enriched library samples may also be qualitatively analyzed using gel electrophoresis. Sample gel electrophoresis results are provided in the [Reference](#) section on [page 105](#).

### Expected Results

Each amplicon in the prepared library contains one target insert surrounded by sequence motifs required for multiplexed sequencing using the Illumina platform. Amplicons include 50 to 500 bp of target DNA insert and 125 bp of sequencing motifs, as shown in [Figure 8](#).



**Figure 8** Content of HaloPlex-enriched target amplicons. Each amplicon contains one target insert (blue) surrounded by the Illumina paired-end sequencing elements (yellow), the sample index (green) and the library bridge PCR primers (gray).

The amplicons should range from 175 to 625 bp in length, and amplicons in the 175 to 625 bp size range should be included for quantitation of the enriched target DNA in each sample. Any spurious DNA products outside of this size range in any sample should be excluded from the target DNA quantitation results.

### 3 Sample Preparation

#### Step 7. Validate enrichment and quantify enriched target DNA

##### Option 1: Analysis using the 2100 Bioanalyzer System

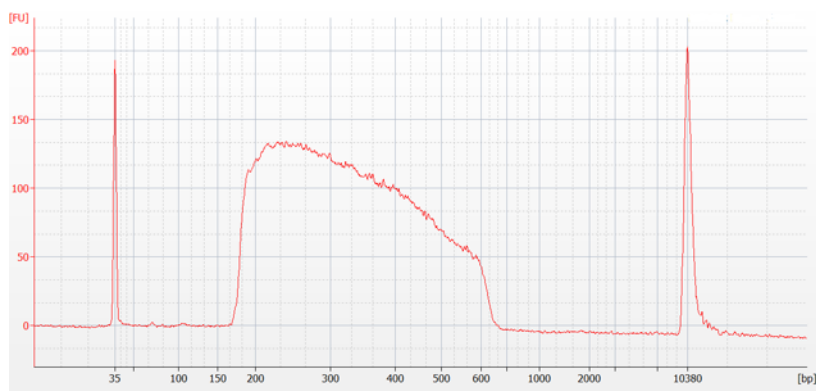
Use a Bioanalyzer High Sensitivity DNA Assay kit and the 2100 Bioanalyzer instrument with 2100 Expert Software (version B.02.07 or higher required to run the High Sensitivity Kit). See the reagent kit guide for general Bioanalyzer instrument and assay setup instructions.

- 1 Prepare the chip, samples and ladder as instructed in the reagent kit guide, using 1  $\mu\text{L}$  of enriched library sample for the analysis.
- 2 Load the prepared chip into the 2100 Bioanalyzer instrument and start the run within five minutes after preparation.
- 3 Analyze the electropherogram for each sample according to the analysis guidelines on [page 84](#).

See [Figure 9](#) for a sample Bioanalyzer electropherogram.

#### NOTE

If the concentration determined by Bioanalyzer analysis is  $> 10 \text{ ng}/\mu\text{L}$ , repeat the analysis using a 1:10 dilution of the sample. Dilute 1  $\mu\text{L}$  of the sample in 9  $\mu\text{L}$  of 10 mM Tris, 1 mM EDTA and then mix well by vortexing at 2000 rpm on the IKA vortex supplied with the Bioanalyzer before analyzing the diluted sample.



**Figure 9** Validation of HaloPlex enrichment by 2100 Bioanalyzer system analysis.

## Step 7. Validate enrichment and quantify enriched target DNA

**Option 2: Analysis using the 2200 TapeStation**

Use a High Sensitivity D1000 ScreenTape (p/n 5067-5584) and reagent kit (p/n 5067-5585) to analyze the enriched library samples. For more information to do this step, see the *2200 TapeStation User Manual*.

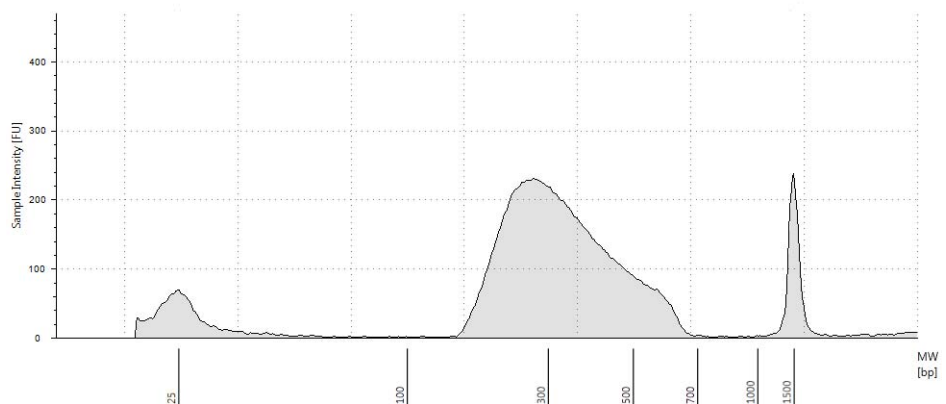
- 1 Prepare the TapeStation samples as instructed in the *2200 TapeStation User Manual*. Use 2  $\mu$ L of each enriched library sample diluted with 2  $\mu$ L of High Sensitivity D1000 sample buffer in separate wells of a tube strip for the analysis.

**CAUTION**

Make sure that you thoroughly mix the combined DNA sample and High Sensitivity D1000 sample buffer on a vortex mixer for 5 seconds for accurate results.

- 2 Load the sample tube strip, the High Sensitivity D1000 ScreenTape, and loading tips into the 2200 TapeStation as instructed in the *2200 TapeStation User Manual*. Start the run.
- 3 Analyze the electropherogram for each sample according to the analysis guidelines on [page 84](#).

See [Figure 10](#) for a sample TapeStation electropherogram.



**Figure 10** Validation of HaloPlex enrichment by 2200 TapeStation analysis.

### 3 Sample Preparation

#### Step 7. Validate enrichment and quantify enriched target DNA

##### **Analysis of Electropherogram Results**

- Check that the electropherogram shows a peak fragment size between approximately 225 to 525 bp.
- Determine the concentration of enriched target DNA in the sample by integration under the peak between 175 and 625 bp. Peaks at <150 bp may be observed, but should be excluded from quantitation.

## Step 8. Pool samples with different indexes for multiplexed sequencing

Use the following guidelines to design your sample pooling strategy:

- Use the Bioanalyzer- or TapeStation-measured concentration of 175-625 bp products in each sample to pool equimolar amounts of differentially indexed samples in order to optimize the use of sequencing capacity.
- The final HaloPlex enrichment pool is ready for direct sequencing using standard Illumina paired-end primers and chemistry on the Illumina HiSeq, MiSeq, or GAIIx platform. See additional guidelines for the MiSeq platform (below) and HiSeq platform ([page 89](#)).
- Use 100 + 100 bp read length paired-end sequencing.
- Sequencing runs must be set up to perform an 8-nt index read. For complete index sequence information, see the [Reference](#) chapter starting on [page 91](#).
- Before aligning reads to the reference genome, trim the reads from Illumina adaptor sequences.

### MiSeq platform sequencing run setup guidelines

#### Setting up a custom Sample Sheet:

- 1 In the IEM software, create a Sample Sheet for the MiSeq platform using the following Workflow selections.
  - Under **Category**, select *Other*.
  - Under **Application**, select *FASTQ Only*.

### 3 Sample Preparation

#### Step 8. Pool samples with different indexes for multiplexed sequencing

- 2 On the **Workflow Parameters** screen, enter the run information, making sure to specify the key parameters highlighted below:

Illumina Experiment Manager

## Sample Sheet Wizard - Workflow Parameters

**FASTQ Only Run Settings**

Reagent Cartridge Barcode\* MSXXXXXXX-300

Sample Prep Kit TruSeq LT

Index Reads ☐ 0 ☒ 1 ☐ 2

Project Name Test Project

Experiment Name Test Experiment

Investigator Name Test

Description Test

Date 2013-01-23

Read Type ☒ Paired End ☐ Single Read

Cycles Read 1 101

Cycles Read 2 101

\* - required field

**FASTQ Only Workflow-Specific Settings**

☐ Custom Primer for Read 1

☐ Custom Primer for Index

☐ Custom Primer for Read 2

☒ Use Adapter Trimming

Cancel Back Next

## Step 8. Pool samples with different indexes for multiplexed sequencing

- 3 Using the **Sample Plate Wizard**, set up a New Plate, entering the required information for each sample to be sequenced. In the **Index 1(17)** column of the **TrueSeq LT Assay Plate** table, assign each sample to any of the Illumina 17 indexes. The index will be corrected to a HaloPlex index at a later stage.

Sample Plate Wizard - Plate Samples

TruSeq LT Assay Plate

Table Plate Plate Graphic ☐ indicates invalid samples

	Sample ID*	Sample Name	Index1 (I7)*	Sample Project	Description
A01	1	Sample1	A001	ProjectX	Tumor
A02	2	Sample2	A002	ProjectX	Normal
A03	3	Sample3	A003	ProjectX	Tumor

- 4 Finish the sample plate setup tasks and save the sample plate file.
- 5 Using the **Sample Sheet Wizard**, select the samples to include in the run and save the Sample Sheet file.

### Step 8. Pool samples with different indexes for multiplexed sequencing

## CAUTION

Kits with indexing primers supplied in a blue plate include indexing primers A01–H12. See [page 95](#) for index nucleotide sequence information.

Kits with indexing primers supplied in a clear plate include indexing primers 1–96. See [page 99](#) through [page 104](#) for index nucleotide sequence information.

- 1 Open the Sample Sheet file in a text editor. For each sample, select the text for the 6-nucleotide index (highlighted below), and replace with the appropriate 8-nucleotide HaloPlex index sequence.

The screenshot shows the ESR software interface. The top menu bar includes File, ESR, Search, View, Encoding, Language, Settings, Macro, Run, Plugins, Window, and ?.

The main window displays the contents of the file `MS000000-300.csv`. The content is as follows:

```
[Header]
IEMFileVersion,4
Investigator Name,Test
Project Name,Test Project
Experiment Name,Test Experiment
Date,9/28/2012
Workflow,GenerateFASTQ
Application,FASTQ Only
Assay,TruSeq LT
Description,Test
Chemistry,Default

[Reads]
151
151

[Settings]

[Data]
Sample_ID,Sample_Name,Sample_Plate,Sample_Well,I7_Index_ID,index,Sample_Project,Description
1,Sample1,testplate,A01,A001,ATCACG,ProjectX,Tumor
2,Sample2,testplate,A02,A002,CAGTGT,ProjectX,Normal
3,Sample3,testplate,A03,A003,TTAGGC,ProjectX,Tumor
4,Sample4,testplate,A04,A004,TGACCA,ProjectX,Normal
5,Sample5,testplate,A05,A005,ACAGTG,ProjectX,Tumor
6,Sample6,testplate,A06,A006,GCCAAAT,ProjectX,Normal
```

- 2 Save the edited Sample Sheet in an appropriate file location for use in the MiSeq platform run.



## Step 8. Pool samples with different indexes for multiplexed sequencing

**HiSeq platform sequencing run setup guidelines**

Set up sequencing runs to perform an 8-nt index read using the *Cycles* settings shown in [Table 35](#). 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 35** HiSeq platform Run Configuration screen Cycle Number settings \*

Run Segment	Cycle Number
Read 1	100
Index 1 (i7)	9
Index 2 (i5)	0
Read 2	100

\* Settings apply to v3.0 SBS chemistry.

**Sequence analysis resources**

Agilent's SureCall data analysis software is available to simplify the sequencing data analysis workflow after HaloPlex target enrichment. To learn more about this resource and download the SureCall software free of charge, visit [www.agilent.com/genomics/surecall](http://www.agilent.com/genomics/surecall).

### **3 Sample Preparation**

#### **Step 8. Pool samples with different indexes for multiplexed sequencing**



## 4 Reference

Reference Information for Kits with Revised Index Configuration (indexing primers in blue plate) [92](#)

Reference Information for Kits with Original Index Configuration (indexing primers in clear plate) [96](#)

Qualitative analysis of enrichment by gel electrophoresis [105](#)

This chapter contains reference information, including component kit contents, index sequences, and optional gel validation instructions.

### CAUTION

This chapter contains two sets of index sequence and kit content information. **Verify that you are referencing the information appropriate for your kit version before you proceed.**

The first section covers kits with reconfigured indexing primers, typically received December, 2014 or later. The reconfigured primers A01–H12 are supplied in a blue plate. See [page 92](#) through [page 95](#) for details.

The second section covers kits with original indexing primer configuration, typically received before December, 2014. The original configuration includes primers 1–96, supplied in a clear plate. See [page 96](#) through [page 104](#) for details.



# Reference Information for Kits with Revised Index Configuration (indexing primers in blue plate)

If your kit includes indexing primers in a clear plate, instead see [page 96](#) for kit content and indexing primer information.

## Kit Contents-Revised Configuration

The HaloPlex Exome Target Enrichment System (revised index configuration) includes the component kits listed below:

**Table 36** HaloPlex Exome Target Enrichment System Kit Contents

Kit Type	HaloPlex Exome, ILM, Box 1 <sup>*</sup>	HaloPlex Magnetic Beads, Box 2
	Store at –20°C	Store at +4°C
HaloPlex Exome, ILM, 96 Reactions, Automated	5190-8064	5190-5386

<sup>\*</sup> See [Table 37](#) for list of included reagents.

The contents of the HaloPlex Exome Target Enrichment System Box 1 (revised index configuration) included with each kit are detailed in the table below:

**Table 37** HaloPlex Exome Target Enrichment System Box 1 Contents

Included Reagents	Format
Hybridization Solution	bottle
Ligation Solution	bottle
Wash Solution	bottle
Capture Solution	bottle
SSC Buffer	bottle
RE Buffer	bottle
BSA Solution	tube with clear cap
DNA Ligase	tube with red cap
Enrichment Control DNA	tube with orange cap
Primer 1	tube with yellow cap
Primer 2	tube with blue cap
HaloPlex Indexing Primer Cassettes	96-well plate with Indexing Primer Cassettes A01 to H12 (blue plate)*
Enzyme Strip 1	8-well strip tube with green label
Enzyme Strip 2	8-well strip tube with red label
HaloPlex Probe 8-well Strip	8-well strip tube with black label

\* See [Table 38](#) for a plate map.

**4 Reference**  
**Kit Contents-Revised Configuration**

**Table 38 Plate map for HaloPlex Indexing Primer Cassettes A01 through H12 provided in blue plate**

	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 HaloPlex Indexes (indexing primers in blue plate)

The nucleotide sequence of the 8-nucleotide index portion of each HaloPlex Indexing Primer Cassette (revised index configuration) is provided in the table below.

**Table 39 HaloPlex Indexes, for indexing primers provided in blue 96-well plate**

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

If your kit includes indexing primers in a blue plate, instead see [page 92](#) for kit content and indexing primer information.

The HaloPlex Exome Target Enrichment System (original index configuration) includes the following component kits:

Kit Type	HaloPlex Exome, ILM, Box 1 *	HaloPlex Magnetic Beads, Box 2
	Store at –20°C	Store at +4°C
<b>HaloPlex Exome, ILM, 96 Reactions, Automated</b>	5190-6291	5190-5386

## HaloPlex Exome Target Enrichment System Automation Protocol



The contents of the HaloPlex Exome Target Enrichment System Box 1 (original index configuration) included with each kit are detailed in the table below:

**Table 41** HaloPlex Exome Target Enrichment System Box 1 Contents

Included Reagents	Format
Hybridization Solution	bottle
Ligation Solution	bottle
Wash Solution	bottle
Capture Solution	bottle
SSC Buffer	bottle
RE Buffer	bottle
BSA Solution	tube with clear cap
DNA Ligase	tube with red cap
Enrichment Control DNA	tube with orange cap
Primer 1	tube with yellow cap
Primer 2	tube with blue cap
HaloPlex Indexing Primer Cassettes	96-well plate with Indexing Primer Cassettes 1-96 (clear plate)*
Enzyme Strip 1	8-well strip tube with green label
Enzyme Strip 2	8-well strip tube with red label
HaloPlex Probe 8-well Strip	8-well strip tube with black label

\* See [Table 42](#) for a plate map.

## 4 Reference

### Kit Contents-Original Configuration

**Table 42 Plate map for HaloPlex Indexing Primer Cassettes 1-96 provided in clear plate**

	1	2	3	4	5	6	7	8	9	10	11	12
<b>A</b>	Index 1	Index 9	Index 17	Index 25	Index 33	Index 41	Index 49	Index 57	Index 65	Index 73	Index 81	Index 89
<b>B</b>	Index 2	Index 10	Index 18	Index 26	Index 34	Index 42	Index 50	Index 58	Index 66	Index 74	Index 82	Index 90
<b>C</b>	Index 3	Index 11	Index 19	Index 27	Index 35	Index 43	Index 51	Index 59	Index 67	Index 75	Index 83	Index 91
<b>D</b>	Index 4	Index 12	Index 20	Index 28	Index 36	Index 44	Index 52	Index 60	Index 68	Index 76	Index 84	Index 92
<b>E</b>	Index 5	Index 13	Index 21	Index 29	Index 37	Index 45	Index 53	Index 61	Index 69	Index 77	Index 85	Index 93
<b>F</b>	Index 6	Index 14	Index 22	Index 30	Index 38	Index 46	Index 54	Index 62	Index 70	Index 78	Index 86	Index 94
<b>G</b>	Index 7	Index 15	Index 23	Index 31	Index 39	Index 47	Index 55	Index 63	Index 71	Index 79	Index 87	Index 95
<b>H</b>	Index 8	Index 16	Index 24	Index 32	Index 40	Index 48	Index 56	Index 64	Index 72	Index 80	Index 88	Index 96

## Nucleotide Sequences of HaloPlex Indexes (indexing primers in clear plate)

The nucleotide sequence of the 8-nucleotide index portion of each HaloPlex Indexing Primer Cassette (original index configuration) is provided in the [Table 43](#) to [Table 48](#) below.

**Table 43** HaloPlex Indexes 1-16

Index Number	Sequence
1	AACGTGAT
2	AAACATCG
3	ATGCCTAA
4	AGTGGTCA
5	ACCACTGT
6	ACATTGGC
7	CAGATCTG
8	CATCAAGT
9	CGCTGATC
10	ACAAGCTA
11	CTGTAGCC
12	AGTACAAG
13	AACAACCA
14	AACCGAGA
15	AACGCTTA
16	AAGACGGA

## 4 Reference

### Nucleotide Sequences of HaloPlex Indexes (indexing primers in clear plate)

**Table 44** HaloPlex Indexes 17-32

Index Number	Sequence
17	AAGGTACA
18	ACACAGAA
19	ACAGCAGA
20	ACCTCCAA
21	ACGCTCGA
22	ACGTATCA
23	ACTATGCA
24	AGAGTCAA
25	AGATCGCA
26	AGCAGGAA
27	AGTCACTA
28	ATCCTGTA
29	ATTGAGGA
30	CAACCACA
31	CAAGACTA
32	CAATGGAA

## Nucleotide Sequences of HaloPlex Indexes (indexing primers in clear plate)

**Table 45** HaloPlex Indexes 33-48

Index Number	Sequence
33	CACTTCGA
34	CAGCGTTA
35	CATACCAA
36	CCAGTTCA
37	CCGAAGTA
38	CCGTGAGA
39	CCTCCTGA
40	CGAACTTA
41	CGACTGGA
42	CGCATACA
43	CTCAATGA
44	CTGAGCCA
45	CTGGCATA
46	GAATCTGA
47	GACTAGTA
48	GAGCTGAA

Nucleotide Sequences of HaloPlex Indexes (indexing primers in clear plate)

**Table 46** HaloPlex Indexes 49-64

Index Number	Sequence
49	GATAGACA
50	GCCACATA
51	GCGAGTAA
52	GCTAACGA
53	GCTCGGTA
54	GGAGAACA
55	GGTGCGAA
56	GTACGCAA
57	GTCGTAGA
58	GTCTGTCA
59	GTGTTCTA
60	TAGGATGA
61	TATCAGCA
62	TCCGTCTA
63	TCTTCACA
64	TGAAGAGA

## Nucleotide Sequences of HaloPlex Indexes (indexing primers in clear plate)

**Table 47** HaloPlex Indexes 65-80

Index Number	Sequence
65	TGGAACAA
66	TGGCTTCA
67	TGGTGGTA
68	TTCACGCA
69	AACTCACC
70	AAGAGATC
71	AAGGACAC
72	AATCCGTC
73	AATGTTGC
74	ACACGACC
75	ACAGATTC
76	AGATGTAC
77	AGCACCTC
78	AGCCATGC
79	AGGCTAAC
80	ATAGCGAC

Nucleotide Sequences of HaloPlex Indexes (indexing primers in clear plate)

**Table 48** HaloPlex Indexes 81-96

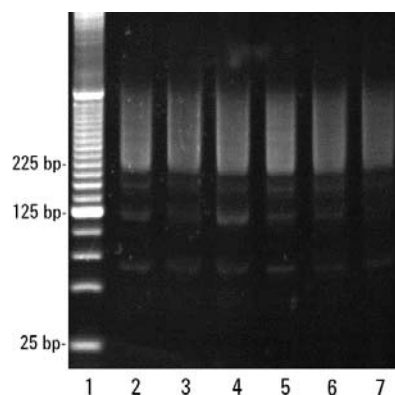
Index Number	Sequence
81	ATCATTCC
82	ATTGGCTC
83	CAAGGAGC
84	CACCTTAC
85	CCATCCTC
86	CCGACAAC
87	CCTAATCC
88	CCTCTATC
89	CGACACAC
90	CGGATTGC
91	CTAAGGTC
92	GAACAGGC
93	GACAGTGC
94	GAGTTAGC
95	GATGAATC
96	GCCAAGAC



## Qualitative analysis of enrichment by gel electrophoresis

Enrichment products may be qualitatively analyzed by gel electrophoresis. Analyze 5  $\mu$ L of each enriched library sample (enriched ECD sample or experimental enriched libraries) by electrophoresis on a Novex 6% polyacrylamide TBE pre-cast gel. See [page 42](#) for additional gel analysis protocol recommendations.

Successful enrichment is indicated by the presence of a smear of amplicons from approximately 225 to 525 bp in each enrichment library lane. For some probe designs, low molecular weight (<150 bp) bands may also be visible, but should not be included in enriched sample quantitation. See [Figure 11](#) for a sample gel analysis image.



**Figure 11** Validation of HaloPlex enrichment process by gel electrophoresis. Lane 1: 25-bp DNA ladder, Lanes 2-7: enriched library samples.

**www.agilent.com**

## **In This Book**

This guide contains information to run the HaloPlex Exome Target Enrichment System automation protocol.

© Agilent Technologies, Inc. 2015

Version C0, July 2015



G9906-90020



**Agilent Technologies**