



# **Preparation of Dual Indexed Libraries using SureSelect<sup>XT</sup> Low Input Target Enrichment System**

**For Illumina Multiplexed  
Sequencing Platforms**

## **Protocol**

Version E0, September 2022

**SureSelect platform manufactured with Agilent  
SurePrint Technology**

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



**Agilent Technologies**

## Notices

© Agilent Technologies, Inc. 2018-2022

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

G9703-90050

### Edition

Version E0, September 2022

Printed in USA

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

### Acknowledgment

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

### Technical Support

#### For US and Canada

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

Or send an e-mail to:

[ngs.support@agilent.com](mailto:ngs.support@agilent.com)

#### For all other regions

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

[www.agilent.com/en/contact-us/page](http://www.agilent.com/en/contact-us/page).

## Warranty

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

## Technology Licenses

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

## Restricted Rights Legend

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

## Notice to Purchaser

This product is provided under an agreement between Bio-Rad Laboratories and Agilent Technologies Inc., and the manufacture, use, sale or import of this product is subject to U.S. Pat. No. 6,670,808, owned by Bio-Rad Laboratories, Inc. Purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in PCR (but not including real-time PCR) in the Research Field (including all Applied Research Fields, including but not limited to forensics, animal testing, and food testing) and in real-time PCR in the Diagnostics and Prognostics Fields. No rights are granted for use of this product for real-time PCR in the Research Field, including all Applied Research Fields (including but not limited to forensics, animal testing and food testing).

## 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 provides an optimized protocol for preparation of dual-indexed, target-enriched Illumina paired-end sequencing libraries using SureSelect<sup>XT</sup> Low Input Reagent Kits and Dual Index P5 Indexed Adaptors 1-96 for ILM.

### 1 Before You Begin

This chapter contains information that you should read and understand before you start an experiment.

### 2 Preparation and Fragmentation of Input DNA

This chapter describes the steps to prepare and fragment gDNA samples, using either mechanical shearing or enzymatic fragmentation, prior to library preparation.

### 3 Library Preparation

This chapter describes the steps to prepare dual-indexed gDNA sequencing libraries for target enrichment.

### 4 Hybridization and Capture

This chapter describes the steps to hybridize and capture the prepared DNA libraries using a SureSelect or ClearSeq Probe.

### 5 Post-Capture Sample Processing for Multiplexed Sequencing

This chapter describes the steps for post-capture amplification and guidelines for sequencing sample preparation.

### 6 Appendix: Using FFPE-derived DNA Samples

This chapter describes the protocol modifications for gDNA isolated from FFPE samples.

### 7 Reference

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

## What's New in Version E0

- Support for SureSelect XT HS Human All Exon V8+UTR and SureSelect XT HS Human All Exon V8+NCV Probes. See [Table 3](#) on page 14 for ordering information.
- Design ID information added to [Table 3](#) on page 14 for pre-designed SureSelect probes.
- Updates to tube cap strip recommendations from domed caps to flat/domed caps based on specific cycler recommendations (see [Table 5](#) on page 16 and related updates on [page 12](#), [page 19](#), and [page 58](#)).
- Update to enzymatic fragmentation instructions in [step 1](#) on [page 29](#) for optional use of 1X Low TE Buffer as solvent for DNA samples.
- Updates to recommended reagent volumes for 24 reaction runs in [Table 15](#) on page 35 and [Table 17](#) on page 36.
- New AmpPure XP bead purification protocol parameter summary tables for experienced users (see [Table 19](#) on page 40, [Table 24](#) on page 45, and [Table 35](#) on page 67).
- Updates to downstream sequencing support information (see [page 78](#) to [page 85](#)) including support for Agilent's AGeNT v3.0 and the included CReaK tool, replacing AGeNT's LocatIt tool (see [page 84](#)).
- Support for use of Agilent's Alissa Reporter software for SureSelect XT HS/Low Input DNA library sequence pre-processing and human germline DNA variant analysis (see [page 78](#) and [page 83](#)).
- Updates to dual index information on [page 96](#) through [page 100](#) to clarify P5 index sequence orientation usage.
- Update to *Troubleshooting* on [page 104](#) on thermal cycler block configuration requirements for efficient heating of SureSelect Wash Buffer 2.
- Update to “[Notice to Purchaser](#)” on page 2.

## What's New in Version D1

- Support for SureSelect XT HS Human All Exon V8 Probe (see [Table 3](#) on page 14)
- Updates to the “[Hybridization and Capture](#)” chapter on [page 53](#) through [page 59](#), including updates to [Table 26](#) on page 55 and additional minor updates throughout the chapter.
- Update to *Note* on [page 26](#) and new footnote to [Table 12](#) on page 30 on impacts of initial FFPE DNA sample fragment size on final library fragment size distribution.

## What's New in Version D0

- Addition of optional Enzymatic DNA Fragmentation protocol (see [page 29](#) to [page 31](#) for the protocol; also see [Table 39](#) on page 88, [Table 43](#) on page 92, and Troubleshooting on [page 101](#))
- New chapter “[Preparation and Fragmentation of Input DNA](#)” starting on [page 21](#) (includes input DNA preparation and mechanical shearing information previously found in “Sample Preparation” chapter)
- Addition of hybridization temperature considerations for probes designed for use with the SureSelect XT system (see footnote to [Table 26](#) on page 55)
- Minor updates to instructions in the “[Hybridization and Capture](#)” chapter, including *Note* on [page 53](#), [step 3](#) on [page 54](#), [step 5](#) on [page 56](#), and [step 1](#) on [page 60](#)
- Updates to downstream sequencing platform and kit support information ([Table 37](#) on page 77, [Table 53](#) on page 99 and [Table 54](#) on page 100)
- Addition of small volume spectrophotometer to [Table 5](#) on page 16
- Update to description of flat strip caps in [Table 7](#) on page 19

# Content

<b>1</b>	<b>Before You Begin</b>	<b>9</b>
	Overview of the Workflow	10
	Procedural Notes	12
	Safety Notes	12
	Materials Required	13
	Optional Materials	19
<b>2</b>	<b>Preparation and Fragmentation of Input DNA</b>	<b>21</b>
	Step 1. Prepare and analyze quality of genomic DNA samples	22
	Preparation of high-quality gDNA from fresh biological samples	22
	Preparation and qualification of gDNA from FFPE samples	23
	Step 2. Fragment the DNA	26
	Method 1: Mechanical DNA Shearing using Covaris	26
	Method 2: Enzymatic DNA Fragmentation	29
<b>3</b>	<b>Library Preparation</b>	<b>33</b>
	Step 1. Repair and dA-Tail the DNA ends	34
	Step 2. Ligate the P5-indexed adaptor	38
	Step 3. Purify the sample using AMPure XP beads	40
	Step 4. Amplify the adaptor-ligated library	42
	Step 5. Purify the amplified library with AMPure XP beads	45
	Step 6. Assess quality and quantity	47

## Contents

<b>4</b>	<b>Hybridization and Capture</b>	<b>53</b>
	Step 1. Hybridize DNA samples to the probe	54
	Step 2. Prepare streptavidin-coated magnetic beads	59
	Step 3. Capture the hybridized DNA using streptavidin-coated beads	60
<b>5</b>	<b>Post-Capture Sample Processing for Multiplexed Sequencing</b>	<b>63</b>
	Step 1. Amplify the captured libraries	64
	Step 2. Purify the amplified captured libraries using AMPure XP beads	67
	Step 3. Assess sequencing library DNA quantity and quality	69
	Step 4. Pool samples for multiplexed sequencing	74
	Step 5. Prepare sequencing samples	76
	Step 6. Do the sequencing run and analyze the data	78
	Sequence analysis resources	83
<b>6</b>	<b>Appendix: Using FFPE-derived DNA Samples</b>	<b>87</b>
	Protocol modifications for FFPE Samples	88
	Methods for FFPE Sample Qualification	88
	Sequencing Output Recommendations for FFPE Samples	89
<b>7</b>	<b>Reference</b>	<b>91</b>
	Kit Contents	92
	Dual Indexing Primer Information	96
	Troubleshooting Guide	101
	Quick Reference Protocol	106





# 1 Before You Begin

Overview of the Workflow	10
Procedural Notes	12
Safety Notes	12
Materials Required	13
Optional Materials	19

Make sure you have the most current protocol. Go to [genomics.agilent.com](http://genomics.agilent.com) and search for G9703-90050.

To prepare dual-indexed libraries for Agilent SureSelect Cancer All-In-One assays, use the protocols detailed in this publication, while implementing the considerations provided in the SureSelect Cancer All-In-One Target Enrichment Product Overview Guide (publication [G9702-90100](#)).

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

## NOTE

This protocol differs from the Illumina Multiplexed Paired-End sequencing manual and other SureSelect protocols at several steps. Make sure to use P5 Indexed Adaptors instead of the Adaptor Oligo Mix and pay close attention to the primers used for each amplification step and the blocking agents used during hybridization.

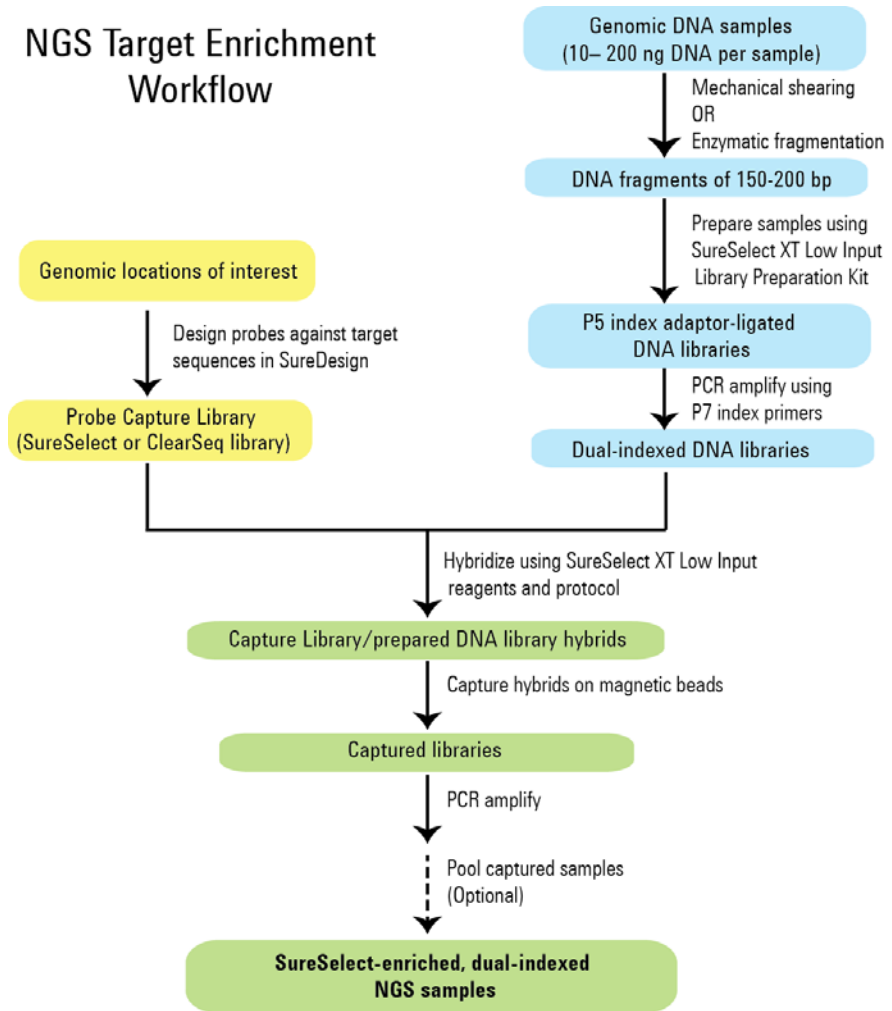
## NOTE

Agilent guarantees performance and provides technical support for the SureSelect reagents required for this workflow only when used as directed in this Protocol.



## Overview of the Workflow

The dual indexing SureSelect<sup>XT</sup> Low Input with Dual Indexing workflow is summarized in Figure 1. The estimated time requirements for each step are summarized in Table 1.



**Figure 1** Overall dual-indexed, target-enriched sequencing sample preparation workflow.

**Table 1** Estimated time requirements (up to 16 sample run size)

Step	Time
Library Preparation	3.5 hours
Hybridization and Capture	3.5 hours
Post-capture amplification	1 hour
QC using Bioanalyzer or TapeStation platform and sample pooling	1.5 hours

## Procedural Notes

- To prevent contamination of reagents by nucleases, always wear powder-free laboratory gloves and use dedicated solutions and pipettors with nuclease-free aerosol-resistant tips.
- Use best-practices to prevent PCR product contamination of samples throughout the workflow:
  - 1 Assign separate pre-PCR and post-PCR work areas and use dedicated equipment, supplies, and reagents in each area. In particular, never use materials designated to post-PCR work areas for pre-PCR segments of the workflow.
  - 2 Maintain clean work areas. Clean pre-PCR surfaces that pose the highest risk of contamination daily using a 10% bleach solution.
  - 3 Always use dedicated pre-PCR pipettors with nuclease-free aerosol-resistant tips to pipette dedicated pre-PCR solutions.
  - 4 Wear powder-free gloves. Use good laboratory hygiene, including changing gloves after contact with any potentially-contaminated surfaces.
- For each protocol step that requires removal of tube cap strips, reseal the tubes with a fresh strip of caps. Reuse of strip caps can cause sample loss, sample contamination, or imprecision in sample temperatures during thermal cycler incubation steps.
- In general, follow Biosafety Level 1 (BSL1) safety rules.
- Possible stopping points, where samples may be stored at  $-20^{\circ}\text{C}$ , are marked in the protocol. Do not subject the samples to multiple freeze/thaw cycles.

## Safety Notes

### CAUTION

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

## Materials Required

Materials required to complete the SureSelect<sup>XT</sup> Low Input Dual Indexing protocol will vary based on the following considerations:

- DNA sample type: high-quality gDNA derived from fresh/fresh-frozen samples vs. FFPE-derived gDNA samples
- DNA fragmentation method used in workflow: mechanical (Covaris-mediated) shearing vs. enzymatic fragmentation

Materials listed in [Table 2](#), plus a compatible target enrichment Probe (selected from [Table 3](#) on page 14) are required for all Dual Indexing workflows. Refer to [Table 4](#) through [Table 6](#) for additional materials needed to complete the protocols according to your DNA sample type/fragmentation method.

**Table 2** SureSelect Reagents for Dual Indexing Workflows

Description	96 Reaction Kit Part Number*
SureSelect XT Low Input Reagent Kit with Index Primers 1–96 for Illumina (ILM) platform†	G9703A
SureSelect XT Low Input Dual Index P5 Indexed Adaptors 1-96 for ILM	5191-4056

\* 96-reaction kits contain enough reagents for 4 runs containing 24 samples per run.

† Compatible with HiSeq, MiSeq, NextSeq 500, and NovaSeq 6000 platforms.

## 1 Before You Begin

### Materials Required

**Table 3** Compatible Probes\*

Probe Capture Library	Design ID	Part Number/Ordering Information (96 Reactions)
<b>Pre-designed Probes</b>		<b>96 Reactions</b>
SureSelect XT HS Human All Exon V8	S33266340	5191-6874
SureSelect XT HS Human All Exon V8+UTR	S33613271	5191-7402
SureSelect XT HS Human All Exon V8+NCV	S33699751	5191-7408
SSel XT HS and XT Low Input Human All Exon V7	S31285117	5191-4029
SureSelect XT Human All Exon V6	S07604514	5190-8864
SureSelect XT Human All Exon V6 + UTRs	S07604624	5190-8882
SureSelect XT Clinical Research Exome V2	S30409818	5190-9492
ClearSeq Comprehensive Cancer XT	0425761	5190-8012
<b>Custom Probes†</b>		
SureSelect Custom Tier1 1–499 kb	Please visit the <a href="#">SureDesign website</a> to design Custom SureSelect probes and obtain ordering information. Contact the SureSelect support team (see <a href="#">page 2</a> ) or your local representative if you need assistance. Custom probes are also available in a 480 Reaction package.	
SureSelect Custom Tier2 0.5–2.9 Mb		
SureSelect Custom Tier3 3–5.9 Mb		
SureSelect Custom Tier4 6–11.9 Mb		
SureSelect Custom Tier5 12–24 Mb		
<b>Pre-designed Probes customized with additional <i>Plus</i> custom content</b>		
SSel XT HS and XT Low Input Human All Exon V7 Plus 1	Please visit the <a href="#">SureDesign website</a> to design the customized <i>Plus</i> content and obtain ordering information. Contact the SureSelect support team (see <a href="#">page 2</a> ) or your local representative if you need assistance.	
SSel XT HS and XT Low Input Human All Exon V7 Plus 2		
SureSelect XT Human All Exon V6 Plus 1		
SureSelect XT Human All Exon V6 Plus 2		
SureSelect XT Clinical Research Exome V2 Plus 1		
SureSelect XT Clinical Research Exome V2 Plus 2		

\* Protocols in this document are also compatible with bundled SureSelect XT Low Input Reagent Kits + Target Enrichment Probes, ordered using p/n G9707A-S. See [page 95](#) for more information.

† Custom Probes designed August 2020 or later are produced using an updated manufacturing process; design-size Tier is shown on labeling for these products. Custom Probes designed and ordered prior to August 2020 may be reordered, with these probes produced using the legacy manufacturing process; design-size Tier is not shown on labeling for the legacy products. Custom Probes of both categories use the same optimized target enrichment protocols detailed in this publication.

**Table 4** Required Reagents--All Sample Types/Fragmentation Methods

Description	Vendor and part number
AMPure XP Kit	Beckman Coulter Genomics
5 ml	p/n A63880
60 ml	p/n A63881
450 ml	p/n A63882
Dynabeads MyOne Streptavidin T1	Thermo Fisher Scientific
2 ml	p/n 65601
10 ml	p/n 65602
50 ml	p/n 65604D
1X Low TE Buffer (10 mM Tris-HCl, pH 7.5-8.0, 0.1 mM EDTA)	Thermo Fisher Scientific p/n 12090-015, or equivalent
100% Ethanol (Ethyl Alcohol, 200 proof)	Millipore p/n EX0276
Qubit BR dsDNA Assay Kit	Thermo Fisher Scientific
100 assays	p/n Q32850
500 assays	p/n Q32853
Nuclease-free Water (not DEPC-treated)	Thermo Fisher Scientific p/n AM9930

## 1 Before You Begin

### Materials Required

#### CAUTION

Sample volumes exceed 0.2 ml in certain steps of this protocol. Make sure that the plasticware used with the selected thermal cycler holds  $\geq 0.25$  ml per well.

**Table 5** Required Equipment--All Sample Types/Fragmentation Methods

Description	Vendor and part number
Thermal Cycler with 96-well, 0.2 ml block	Various suppliers
Plasticware compatible with the selected thermal cycler: 96-well plates or 8-well strip tubes Tube cap strips (flat or domed, based on cycler/lid requirements)*	Consult the thermal cycler manufacturer's recommendations
Qubit Fluorometer	Thermo Fisher Scientific p/n Q33238
Qubit Assay Tubes	Thermo Fisher Scientific p/n Q32856
DNA LoBind Tubes, 1.5-ml PCR clean, 250 pieces	Eppendorf p/n 022431021 or equivalent
Microcentrifuge	Eppendorf microcentrifuge, model 5417C or equivalent
Plate or strip tube centrifuge	Labnet International MPS1000 Mini Plate Spinner, p/n C1000 (requires adapter, p/n C1000-ADAPT, for use with strip tubes) or equivalent
96-well plate mixer	Eppendorf ThermoMixer C, p/n 5382000023 and Eppendorf SmartBlock 96 PCR, p/n 5306000006, or equivalent
Small-volume spectrophotometer	NanoDrop 2000, Thermo Fisher Scientific p/n ND-2000 or equivalent
Multichannel pipette	Rainin Pipet-Lite Multi Pipette or equivalent
Single channel pipettes (10-, 20-, 200-, and 1000- $\mu$ l capacity)	Rainin Pipet-Lite Pipettes or equivalent
Sterile, nuclease-free aerosol barrier pipette tips	general laboratory supplier
Vortex mixer	general laboratory supplier



**Table 5** Required Equipment--All Sample Types/Fragmentation Methods

Description	Vendor and part number
DNA Analysis Platform and Consumables <sup>†</sup>	
Agilent 2100 Bioanalyzer Instrument	Agilent p/n G2939BA
Agilent 2100 Expert SW Laptop Bundle (optional)	Agilent p/n G2953CA
DNA 1000 Kit	Agilent p/n 5067-1504
High Sensitivity DNA Kit	Agilent p/n 5067-4626
OR	
Agilent 4200 TapeStation	Agilent p/n G2991AA/G2992AA
96-well sample plates	Agilent p/n 5042-8502
96-well plate foil seals	Agilent p/n 5067-5154
8-well tube strips	Agilent p/n 401428
8-well tube strip caps	Agilent p/n 401425
D1000 ScreenTape	Agilent p/n 5067-5582
D1000 Reagents	Agilent p/n 5067-5583
High Sensitivity D1000 ScreenTape	Agilent p/n 5067-5584
High Sensitivity D1000 Reagents	Agilent p/n 5067-5585
Magnetic separator	Thermo Fisher Scientific p/n 12331D or equivalent <sup>‡</sup>
Ice bucket	general laboratory supplier
Powder-free gloves	general laboratory supplier

\* Consult the thermal cycler manufacturer's recommendations for use of either flat or domed strip caps and for any accessories (e.g., compression mats) required for optimal performance with the selected caps. Ensure that the combination of selected thermal cycler and plasticware provides complete sealing of sample wells and optimal contact between the instrument heated lid and vial cap for heat transfer.

† DNA samples may also be analyzed using the Agilent 5200 Fragment Analyzer, p/n M5310AA, and associated NGS Fragment Kits (DNF-473-0500 and DNF-474-0500). Implement any sample dilution instructions provided in protocols in this document, and then follow the assay instructions provided for each NGS Fragment Kit.

‡ Select a magnetic separator configured to collect magnetic particles on one side of each well. Do not use a magnetic separator configured to collect the particles in a ring formation.

## 1 Before You Begin

### Materials Required

**Table 6** Additional Required Materials based on DNA Sample Type/Fragmentation Method

Description	Vendor and Part Number
<b>Required for preparation of high-quality DNA samples (not required for FFPE DNA sample preparation)</b>	
High-quality gDNA purification system, for example:	
QIAamp DNA Mini Kit	Qiagen
50 Samples	p/n 51304
250 Samples	p/n 51306
<b>Required for preparation of FFPE DNA samples (not required for high-quality DNA sample preparation)</b>	
QIAamp DNA FFPE Tissue Kit, 50 Samples	Qiagen p/n 56404
Deparaffinization Solution	Qiagen p/n 19093
FFPE DNA integrity assessment system:	
Agilent NGS FFPE QC Kit	Agilent
16 reactions	p/n G9700A
96 reactions	p/n G9700B
<b>OR</b>	
TapeStation Genomic DNA Analysis Consumables:	Agilent
Genomic DNA ScreenTape	p/n 5067-5365
Genomic DNA Reagents	p/n 5067-5366
<b>Required for mechanical shearing of DNA samples (not required for workflows with enzymatic fragmentation)</b>	
Covaris Sample Preparation System	Covaris model E220
Covaris microTUBE sample holders	Covaris p/n 520045
<b>Required for enzymatic fragmentation of DNA samples (not required for workflows with mechanical shearing)</b>	
SureSelect Enzymatic Fragmentation Kit	Agilent
	p/n 5191-4079 (16 reactions)
	p/n 5191-4080 (96 reactions)

## Optional Materials

**Table 7** Supplier Information for Optional Materials

Description	Vendor and Part Number	Purpose
Tween 20	Sigma-Aldrich p/n P9416-50ML	Sequencing library storage (see <a href="#">page 73</a> )
MicroAmp Clear Adhesive Film	Thermo Fisher Scientific p/n 4311971	Improved sealing for flat strip caps
PlateLoc Thermal Microplate Sealer with Small Hotplate and Peelable Aluminum Seal for PlateLoc Sealer	Please contact the SureSelect support team (see <a href="#">page 2</a> ) or your local representative for ordering information	Sealing wells for protocol steps performed inside or outside of the thermal cycler

**1 Before You Begin**  
Optional Materials



## 2 Preparation and Fragmentation of Input DNA

Step 1. Prepare and analyze quality of genomic DNA samples	22
Preparation of high-quality gDNA from fresh biological samples	22
Preparation and qualification of gDNA from FFPE samples	23
Step 2. Fragment the DNA	26
Method 1: Mechanical DNA Shearing using Covaris	26
Method 2: Enzymatic DNA Fragmentation	29

This chapter describes the steps to prepare, quantify, qualify, and fragment input DNA samples prior to SureSelect XT Low Input library preparation and target enrichment. Protocols are provided for two alternative methods of DNA fragmentation—mechanical shearing or enzymatic DNA fragmentation.

The library preparation protocol is compatible with both high-quality gDNA prepared from fresh or fresh-frozen samples and lower-quality DNA prepared from FFPE samples. Modifications required for FFPE samples are included throughout the protocol steps. For a summary of modifications for FFPE samples see [Chapter 6](#), “Appendix: Using FFPE-derived DNA Samples” on [page 87](#).

The protocol requires 10 ng to 200 ng of input DNA, with adjustments to DNA input amount or quantification method required for some FFPE samples. For optimal sequencing results, use the maximum amount of input DNA available within the recommended range.



## 2 Preparation and Fragmentation of Input DNA

### Step 1. Prepare and analyze quality of genomic DNA samples

## Step 1. Prepare and analyze quality of genomic DNA samples

#### NOTE

If you are preparing DNA samples for an Agilent SureSelect Cancer All-In-One assay, use the following modifications to the gDNA sample preparation instructions in this section:

- Where required for your experimental design, make sure to prepare reference DNA sample(s) alongside your experimental samples
- Use at least 50 ng input gDNA for best results

See publication [G9702-90100](#) for more information.

---

### Preparation of high-quality gDNA from fresh biological samples

- 1 Prepare high-quality gDNA using a suitable purification system, such as Qiagen's QIAamp DNA Mini Kit, following the manufacturer's protocol. The protocol requires 10 ng to 200 ng DNA input.

#### NOTE

Make sure genomic DNA samples are of high quality with an OD 260/280 ratio ranging from 1.8 to 2.0.

---

- 2 Use the Qubit BR dsDNA Assay Kit to determine the concentration of each gDNA sample. Follow the manufacturer's instructions for the instrument and assay kit.

Additional qualification of DNA samples is not required for DNA derived from fresh biological samples. Proceed to “[Step 2. Fragment the DNA](#)” on page 26.

## Preparation and qualification of gDNA from FFPE samples

- 1 Prepare gDNA from FFPE tissue sections using Qiagen's QIAamp DNA FFPE Tissue Kit and Qiagen's Deparaffinization Solution, following the manufacturer's protocol. Elute the final gDNA samples from the MinElute column in two rounds, using 30  $\mu$ l Buffer ATE in each round, for a final elution volume of approximately 60  $\mu$ l.

### NOTE

If tissue lysis appears incomplete after one hour of digestion with Proteinase K, add an additional 10  $\mu$ l of Proteinase K and continue incubating at 56°C, with periodic mixing, for up to three hours.

Store the gDNA samples on ice for same-day library preparation, or at -20°C for later processing.

- 2 Assess the quality (DNA integrity) for each FFPE DNA sample using one of the methods below.

### Option 1: Qualification using the Agilent NGS FFPE QC Kit (Recommended Method)

The Agilent NGS FFPE QC Kit provides a qPCR-based assay for DNA sample integrity determination. Results include a  $\Delta\Delta$ Cq DNA integrity score and the precise quantity of amplifiable DNA in the sample, allowing direct normalization of DNA input for each sample. DNA input recommendations based on  $\Delta\Delta$ Cq scores for individual samples are summarized in [Table 8](#).

- a Use the Qubit BR dsDNA Assay Kit to determine the concentration of each gDNA sample. Follow the manufacturer's instructions for the instrument and assay kit.
- b Remove a 1  $\mu$ l aliquot of the FFPE gDNA sample for analysis using the Agilent NGS FFPE QC Kit to determine the  $\Delta\Delta$ Cq DNA integrity score. See the kit user manual (G9700-90000) at [www.agilent.com](http://www.agilent.com) for more information.
- c For all samples with  $\Delta\Delta$ Cq DNA integrity score  $\leq 1$ , use the Qubit-based gDNA concentration determined in [step a](#), above, to determine volume of input DNA needed for the protocol.

## 2 Preparation and Fragmentation of Input DNA

### Preparation and qualification of gDNA from FFPE samples

- d For all samples with  $\Delta\Delta Cq$  DNA integrity score  $>1$ , use the qPCR-based concentration of amplifiable gDNA, reported by the Agilent NGS FFPE QC Kit results, to determine amounts of input DNA for the protocol.

**Table 8** SureSelect XT Low Input DNA input modifications based on  $\Delta\Delta Cq$  DNA integrity score

Protocol Parameter	non-FFPE Samples	FFPE Samples	
		$\Delta\Delta Cq \leq 1^*$	$\Delta\Delta Cq > 1$
DNA input for Library Preparation	10 ng to 200 ng DNA, based on Qubit Assay	10 ng to 200 ng DNA, based on Qubit Assay	10 ng to 200 ng of amplifiable DNA, based on qPCR quantification

\* FFPE samples with  $\Delta\Delta Cq$  scores  $\leq 1$  should be treated like non-FFPE samples for DNA input amount determinations. For samples of this type, make sure to use the DNA concentration determined by the Qubit Assay, instead of the concentration determined by qPCR, to calculate the volume required for 10–200 ng DNA.

#### Option 2: Qualification using Agilent’s Genomic DNA ScreenTape assay DIN score

Agilent’s Genomic DNA ScreenTape assay, used in conjunction with Agilent’s TapeStation, provides a quantitative electrophoretic assay for DNA sample integrity determination. This assay reports a DNA Integrity Number (DIN) score for each sample which is used to estimate the appropriate normalization of DNA input required for low-integrity DNA samples.

- a Use the Qubit BR dsDNA Assay Kit to determine the concentration of each gDNA sample. Follow the manufacturer’s instructions for the instrument and assay kit.
- b Remove a 1  $\mu$ l aliquot of the FFPE gDNA sample and analyze using the Genomic DNA ScreenTape assay. See the user manual at [www.agilent.com](http://www.agilent.com) for more information.
- c Using the DIN score reported for each sample in the Genomic DNA ScreenTape assay, consult [Table 9](#) to determine the recommended amount of input DNA for the sample.



**Table 9** SureSelect XT Low Input DNA input modifications based on DNA Integrity Number (DIN) score

Protocol Parameter	non-FFPE Samples	FFPE Samples		
		DIN > 8*	DIN 3–8	DIN < 3
DNA input for Library Preparation	10 ng to 200 ng DNA, quantified by Qubit Assay	10 ng to 200 ng DNA, quantified by Qubit Assay	Use at least 15 ng for more intact samples and at least 40 ng for less intact samples. Use the maximum amount of DNA available, up to 200 ng, for all samples. Quantify by Qubit Assay.	Use at least 50 ng for more intact samples and at least 100 ng for the least intact samples. Use the maximum amount of DNA available, up to 200 ng, for all samples. Quantify by Qubit Assay.

\* FFPE samples with DIN > 8 should be treated like non-FFPE samples for DNA input amount determinations.

## Step 2. Fragment the DNA

### Method 1: Mechanical DNA Shearing using Covaris

In this step, 50- $\mu$ l gDNA samples are sheared using conditions optimized for either high-quality or FFPE DNA. The target DNA fragment size is 150 to 200 bp.

#### NOTE

This protocol has been optimized using a Covaris model E220 instrument and the 130- $\mu$ l Covaris microTUBE for a target DNA fragment size of 150 to 200 bp. If you wish to use a different Covaris instrument model/sample holder or if your NGS workflow requires a different DNA fragment size (e.g., for translocation detection with the SureSelect Cancer All-In-One assay), consult the manufacturer's recommendations for shearing conditions for the recommended DNA fragment size.

For FFPE DNA samples, initial DNA fragment size may impact the post-shear fragment size distribution, resulting in fragment sizes shorter than the target ranges listed. All FFPE samples should be sheared for 240 seconds (see [Table 10](#) on page 27) to generate fragment ends suitable for library construction. Libraries prepared from FFPE samples should be analyzed using an NGS read length suitable for the final library fragment size distribution.

- 1 Set up the Covaris E220 instrument. Refer to the Covaris instrument user guide for details.
  - a Check that the water in the Covaris tank is filled with fresh deionized water to the appropriate fill line level according to the manufacturer's recommendations for the specific instrument model and sample tube or plate in use.
  - b Check that the water covers the visible glass part of the tube.
  - c On the instrument control panel, push the Degas button. Degas the instrument according to the manufacturer's recommendations, typically 30–60 minutes.
  - d Set the chiller temperature to between 2°C to 5°C to ensure that the temperature reading in the water bath displays 5°C. Consult the manufacturer's recommendations for addition of coolant fluids to prevent freezing.

- 2 Prepare the DNA samples for the run by diluting 10–200 ng of each gDNA sample with 1X Low TE Buffer (10 mM Tris-HCl, pH 7.5-8.0, 0.1 mM EDTA) to a final volume of 50 µl. Vortex well to mix, then spin briefly to collect the liquid. Keep the samples on ice.

**NOTE**

**Do not dilute samples to be sheared using water.** Shearing samples in water reduces the overall library preparation yield and complexity.

- 3 Complete the DNA shearing steps below for each of the gDNA samples.
  - a Transfer the 50-µl DNA sample into a Covaris microTUBE, using a tapered pipette tip to slowly transfer the sample through the pre-split septum of the cap.
  - b Spin the microTUBE for 30 seconds to collect the liquid and to remove any bubbles from the bottom of the tube.
  - c Secure the microTUBE in the tube holder and shear the DNA with the settings in [Table 10](#).

**Table 10** Shear settings for Covaris E-series instrument (SonoLab software v7 or later)

Setting	High-quality DNA	FFPE DNA
Duty Factor	10%	10%
Peak Incident Power (PIP)	175	175
Cycles per Burst	200	200
Treatment Time	2 × 120 seconds	240 seconds
Bath Temperature	2° to 8° C	2° to 8° C

Use the steps below for two-round shearing of **high-quality DNA samples only**:

- Shear for 120 seconds
- Spin the microTUBE for 10 seconds
- Vortex the microTUBE at high speed for 5 seconds
- Spin the microTUBE for 10 seconds
- Shear for additional 120 seconds
- Spin the microTUBE for 10 seconds
- Vortex the microTUBE at high speed for 5 seconds
- Spin the microTUBE for 10 seconds

## 2 Preparation and Fragmentation of Input DNA

### Method 1: Mechanical DNA Shearing using Covaris

- d** After completing the shearing step(s), put the Covaris microTUBE back into the loading and unloading station.
- e** While keeping the snap-cap on, insert a pipette tip through the pre-split septum, then slowly remove the sheared DNA.
- f** Transfer the sheared DNA sample (approximately 50  $\mu$ l) to a 96-well plate or strip tube sample well. Keep the samples on ice.
- g** After transferring the DNA sample, spin the microTUBE briefly to collect any residual sample volume. Transfer any additional collected liquid to the sample well used in [step f](#).

#### NOTE

It is important to avoid loss of input DNA at this step, especially for low-abundance DNA samples. Visually inspect the microTUBE to ensure that all of the sample has been transferred. If droplets remain in the microTUBE, repeat [step g](#).

---

The 50- $\mu$ l sheared DNA samples are now ready for NGS sequencing library preparation, beginning with end repair/dA-tailing. Proceed to “[Library Preparation](#)” on page 33.

#### NOTE

This is not a stopping point in the workflow, and analysis of the sheared samples is not required before they are used for library preparation. Proceed directly to end-repair and dA-tailing.

---

## Method 2: Enzymatic DNA Fragmentation

In this step, gDNA samples are fragmented using Agilent’s SureSelect Enzymatic Fragmentation Kit.

- 1 In wells of a thermal cycler-compatible strip tube or PCR plate, dilute 10 ng to 200 ng of each gDNA sample with nuclease-free water or 1X Low TE Buffer to a final volume of 7  $\mu$ l.

If the DNA concentration is too low to supply the 10–200 ng input amount required for your workflow in 7  $\mu$ l, sample volume may be reduced using a suitable concentration method. Alternatively, see *Troubleshooting* on [page 101](#) for protocol modifications for dilute samples.

- 2 Thaw the vial of 5X SureSelect Fragmentation Buffer on ice, vortex, then keep on ice.
- 3 Preprogram a thermal cycler (with the heated lid ON) with the program in [Table 11](#). Immediately pause the program, and keep paused until samples are loaded in [step 7](#).

**Table 11** Thermal cycler program for enzymatic fragmentation\*

Step	Temperature	Time
Step 1	37°C	Varies—see <a href="#">Table 12</a>
Step 2	65°C	5 minutes
Step 3	4°C	Hold

\* Use a reaction volume setting of 10  $\mu$ l, if required for thermal cycler set up.

Optimal fragmentation conditions may vary based on the NGS read length to be used in the workflow. Refer to [Table 12](#) for the duration at 37°C appropriate for your sample type and required NGS read length.

## 2 Preparation and Fragmentation of Input DNA

### Method 2: Enzymatic DNA Fragmentation

**Table 12** Fragmentation duration based on sample type and NGS read length

NGS read length requirement	Target fragment size	Duration of 37°C incubation step (Table 11)	
		High-quality DNA samples	FFPE DNA samples*
2 × 100 reads	150 to 200 bp	15 minutes	15 minutes
2 × 150 reads	180 to 250 bp	10 minutes	15 minutes

\* For FFPE DNA samples, initial DNA fragment size may impact the post-fragmentation size distribution, resulting in fragment sizes shorter than the target ranges listed in this table. All FFPE samples should be incubated at 37°C for 15 minutes to generate fragment ends suitable for library construction. Libraries prepared from FFPE samples should be analyzed using an NGS read length suitable for the final library fragment size distribution.

- 4 Prepare the appropriate volume of Fragmentation master mix by combining the reagents in Table 13.

Mix well by pipetting up and down 20 times or seal the tube and vortex at high speed for 5–10 seconds. Spin briefly to remove any bubbles and keep on ice.

**Table 13** Preparation of Fragmentation master mix

Reagent	Volume for 1 reaction	Volume for 8 reactions (includes excess)	Volume for 24 reactions (includes excess)
5X SureSelect Fragmentation Buffer (blue cap)	2 µl	18 µl	50 µl
SureSelect Fragmentation Enzyme (green cap)	1 µl	9 µl	25 µl
Total	3 µl	27 µl	75 µl

- 5 Add 3 µl of the Fragmentation master mix to each sample well containing 7 µl of input DNA.
- 6 Mix well by pipetting up and down 20 times or cap the wells and vortex at high speed for 5–10 seconds. Spin the samples briefly.
- 7 Immediately place the plate or strip tube in the thermal cycler and resume the thermal cycling program in Table 11.

- 8 When the program reaches the 4°C Hold step, remove the samples from the thermal cycler, add 40 µl of nuclease-free water to each sample, and place the samples on ice.

The 50-µl reactions are now ready for NGS sequencing library preparation, beginning with end repair/dA-tailing. Proceed to “[Library Preparation](#)” on page 33.

**NOTE**

This is not a stopping point in the workflow, and analysis of the enzymatically-fragmented samples is not required before they are used for library preparation. Proceed directly to end-repair and dA-tailing.

---

## **2 Preparation and Fragmentation of Input DNA**

### **Method 2: Enzymatic DNA Fragmentation**





## 3 Library Preparation

- Step 1. Repair and dA-Tail the DNA ends 34
- Step 2. Ligate the P5-indexed adaptor 38
- Step 3. Purify the sample using AMPure XP beads 40
- Step 4. Amplify the adaptor-ligated library 42
- Step 5. Purify the amplified library with AMPure XP beads 45
- Step 6. Assess quality and quantity 47

The sample preparation protocol is used to prepare DNA libraries for sequencing using the Illumina paired-read platform. For each sample to be sequenced, an individual dual-indexed library is prepared. For an overview of the SureSelect<sup>XT</sup> Low Input target enrichment workflow, see [Figure 1 on page 10](#).

The NGS library preparation protocol that begins here is used for fragmented DNA samples produced by mechanical shearing (as detailed on [page 26 to page 28](#)) or produced by enzymatic fragmentation (as detailed on [page 29 to page 31](#)). Samples produced by either method should contain 10–200 ng of DNA fragments in a volume of 50 µl.



### 3 Library Preparation

#### Step 1. Repair and dA-Tail the DNA ends

## Step 1. Repair and dA-Tail the DNA ends

Protocol steps in this section use the components listed in [Table 14](#). Thaw and mix each component as directed in [Table 14](#) before use.

Remove the AMPure XP beads from cold storage and equilibrate to room temperature in preparation for use on [page 40](#). *Do not freeze the beads at any time.*

**Table 14** Reagents thawed before use in protocol

Kit Component	Storage Location	Thawing Conditions	Mixing Method	Where Used
End Repair-A Tailing Buffer (bottle)	SureSelect XT HS and XT Low Input Library Preparation Kit for ILM (Pre PCR), -20°C	Thaw on ice (may require >20 minutes) then keep on ice	Vortexing	<a href="#">page 36</a>
Ligation Buffer (bottle)	SureSelect XT HS and XT Low Input Library Preparation Kit for ILM (Pre PCR), -20°C	Thaw on ice (may require >20 minutes) then keep on ice	Vortexing	<a href="#">page 35</a>
End Repair-A Tailing Enzyme Mix (orange cap)	SureSelect XT HS and XT Low Input Library Preparation Kit for ILM (Pre PCR), -20°C	Place on ice just before use	Inversion	<a href="#">page 36</a>
T4 DNA Ligase (blue cap)	SureSelect XT HS and XT Low Input Library Preparation Kit for ILM (Pre PCR), -20°C	Place on ice just before use	Inversion	<a href="#">page 35</a>
P5 Indexed Adaptors (green plate)	SureSelect XT Low Input Dual Index P5 Indexed Adaptors 1-96 for ILM, -20°C	Thaw on ice then keep on ice	Vortexing	<a href="#">page 39</a>

To process multiple samples, prepare reagent mixtures with overage at each step, without the DNA sample. Mixtures for preparation of 8 samples and 24 samples (including excess) are shown in each table as examples.

- 1 Before starting the end-repair protocol, prepare the Ligation master mix to allow equilibration to room temperature before use.
  - a Vortex the thawed vial of Ligation Buffer for 15 seconds at high speed to ensure homogeneity.

**CAUTION**

The Ligation Buffer used in this step is viscous. Mix thoroughly by vortexing at high speed for 15 seconds before removing an aliquot for use. When combining with other reagents, mix well by pipetting up and down 15–20 times using a pipette set to at least 80% of the mixture volume or by vortexing at high speed for 10–20 seconds.

Use a flat-top vortex mixer when vortexing strip tubes or plates throughout the protocol. When reagents are mixed by vortexing, visually verify that adequate mixing is occurring.

- b Prepare the appropriate volume of Ligation master mix by combining the reagents in [Table 15](#).

Slowly pipette the Ligation Buffer into a 1.5-ml Eppendorf tube, ensuring that the full volume is dispensed. Slowly add the T4 DNA Ligase, rinsing the enzyme tip with buffer solution after addition. Mix well by slowly pipetting up and down 15–20 times or seal the tube and vortex at high speed for 10–20 seconds. Spin briefly to collect the liquid.

**Keep at room temperature for 30–45 minutes** before use on [page 38](#).

**Table 15** Preparation of Ligation master mix

Reagent	Volume for 1 reaction	Volume for 8 reactions* (includes excess)	Volume for 24 reactions* (includes excess)
Ligation Buffer (bottle)	23 $\mu$ l	207 $\mu$ l	598 $\mu$ l
T4 DNA Ligase (blue cap)	2 $\mu$ l	18 $\mu$ l	52 $\mu$ l
<b>Total</b>	<b>25 <math>\mu</math>l</b>	<b>225 <math>\mu</math>l</b>	<b>650 <math>\mu</math>l</b>

\* The minimum supported run size for 96-reaction kits is 24 samples per run, with kits containing enough reagents for 4 runs of 24 samples each.

### 3 Library Preparation

#### Step 1. Repair and dA-Tail the DNA ends

- 2 Preprogram a thermal cycler (with the heated lid ON) with the program in [Table 16](#) for the End Repair and dA-Tailing steps. Immediately pause the program, and keep paused until samples are loaded in [step 6](#).

**Table 16** Thermal cycler program for End Repair/dA-Tailing \*

Step	Temperature	Time
Step 1	20°C	15 minutes
Step 2	72°C	15 minutes
Step 3	4°C	Hold

\* When setting up the thermal cycling program, use a reaction volume setting of 70  $\mu$ L.

- 3 Vortex the thawed vial of End Repair-A Tailing Buffer for 15 seconds at high speed to ensure homogeneity. Visually inspect the solution; if any solids are observed, continue vortexing until all solids are dissolved.

#### CAUTION

The End Repair-A Tailing Buffer used in this step must be mixed thoroughly by vortexing at high speed for 15 seconds before removing an aliquot for use. When combining with other reagents, mix well either by pipetting up and down 15–20 times using a pipette set to at least 80% of the mixture volume or by vortexing at high speed for 5–10 seconds.

- 4 Prepare the appropriate volume of End Repair/dA-Tailing master mix by combining the reagents in [Table 17](#).

Slowly pipette the End Repair-A Tailing Buffer into a 1.5-ml Eppendorf tube, ensuring that the full volume is dispensed. Slowly add the End Repair-A Tailing Enzyme Mix, rinsing the enzyme tip with buffer solution after addition. Mix well by pipetting up and down 15–20 times or seal the tube and vortex at high speed for 5–10 seconds. Spin briefly to collect the liquid and keep on ice.

**Table 17** Preparation of End Repair/dA-Tailing master mix

Reagent	Volume for 1 reaction	Volume for 8 reactions (includes excess)	Volume for 24 reactions (includes excess)
End Repair-A Tailing Buffer (bottle)	16 $\mu$ l	144 $\mu$ l	416 $\mu$ l
End Repair-A Tailing Enzyme Mix (orange cap)	4 $\mu$ l	36 $\mu$ l	104 $\mu$ l
<b>Total</b>	<b>20 <math>\mu</math>l</b>	<b>180 <math>\mu</math>l</b>	<b>520 <math>\mu</math>l</b>

## Step 1. Repair and dA-Tail the DNA ends

- 5 Add 20  $\mu$ l of the End Repair/dA-Tailing master mix to each sample well containing approximately 50  $\mu$ l fragmented DNA. Mix by pipetting up and down 15–20 times using a pipette set to 60  $\mu$ l or cap the wells and vortex at high speed for 5–10 seconds.
- 6 Briefly spin the samples, then immediately place the plate or strip tube in the thermal cycler and resume the thermal cycling program in [Table 16](#).

### 3 Library Preparation

#### Step 2. Ligate the P5-indexed adaptor

## Step 2. Ligate the P5-indexed adaptor

Before completing this step, assign a dual index pair to each sample. Use indexing pairs made up of the SureSelect XT Low Input Dual Index P5 Indexed Adaptor (green plate) and the SureSelect XT Low Input Index Primer (providing the P7 index; yellow plate) from the same well position of each index plate. Use a different pair of indexes for each sample to be sequenced in the same lane.

The P5 Indexed Adaptors are ligated to DNA libraries in [step 4](#) below. (See [Table 53](#) on page 99 or [Table 54](#) on page 100 for sequences of the 8-bp P5 index portion of the adaptors.) The P7 indexes are introduced during PCR amplification on [page 44](#). (See [Table 52](#) on page 98 for sequences of the 8-bp P7 index portion of the primers.)

- 1 Once the thermal cycler program for End Repair/dA-Tailing reaches the 4°C Hold step, transfer the samples to ice while setting up this step.
- 2 Preprogram a thermal cycler (with the heated lid ON) for the Ligation step with the program in [Table 18](#). Immediately pause the program, and keep paused until samples are loaded in [step 5](#).

**Table 18** Thermal cycler program for Ligation\*

Step	Temperature	Time
Step 1	20°C	30 minutes
Step 2	4°C	Hold

\* Use a reaction volume setting of 100 µl, if required for thermal cycler set up.

- 3 To each end-repaired/dA-tailed DNA sample (approximately 70 µl), add 25 µl of the Ligation master mix that was prepared on [page 35](#) and kept at room temperature. Mix by pipetting up and down at least 10 times using a pipette set to 85 µl or cap the wells and vortex at high speed for 5–10 seconds. Briefly spin the samples.

- 4 Add 5  $\mu$ l of the appropriate SureSelect XT Low Input Dual Index P5 Indexed Adaptor (green plate) to each sample. Use a different adaptor for each sample. Mix by pipetting up and down 15–20 times using a pipette set to 85  $\mu$ l or cap the wells and vortex at high speed for 5–10 seconds.

**NOTE**

Make sure to add the Ligation master mix and the P5 Indexed Adaptor to the samples in separate addition steps as directed in [step 3](#) and [step 4](#) above, mixing after each addition.

- 5 Briefly spin the samples, then immediately place the plate or strip tube in the thermal cycler and resume the thermal cycling program in [Table 18](#).

**Stopping Point** If you do not continue to the next step, seal the sample wells and store overnight at either 4°C or –20°C.

### 3 Library Preparation

#### Step 3. Purify the sample using AMPure XP beads

## Step 3. Purify the sample using AMPure XP beads

In this step, the DNA libraries are purified using AmpPure XP Beads. Critical purification protocol parameters are summarized for experienced users in [Table 19](#).

**Table 19** AMPure XP bead cleanup parameters after adaptor ligation

Parameter	Value
Volume of RT AMPure XP bead suspension added to each sample well	80 $\mu$ l
Final elution solvent and volume	35 $\mu$ l nuclease-free water
Amount of eluted sample transferred to fresh well	Approximately 34.5 $\mu$ l

- 1 Verify that the AMPure XP beads were held at room temperature for at least 30 minutes before use.
- 2 Prepare 400  $\mu$ l of 70% ethanol per sample, plus excess, for use in [step 8](#).

#### NOTE

The freshly-prepared 70% ethanol may be used for subsequent purification steps run on the same day. The complete Library Preparation protocol requires 0.8 ml of fresh 70% ethanol per sample.

- 3 Mix the AMPure XP bead suspension well so that the reagent appears homogeneous and consistent in color.
- 4 Add 80  $\mu$ l of homogeneous AMPure XP beads to each DNA sample (approximately 100  $\mu$ l) in the PCR plate or strip tube. Pipette up and down 15–20 times or cap the wells and vortex at high speed for 5–10 seconds to mix.
- 5 Incubate samples for 5 minutes at room temperature.
- 6 Put the plate or strip tube into a magnetic separation device. Wait for the solution to clear (approximately 5 to 10 minutes).
- 7 Keep the plate or strip tube in the magnetic stand. Carefully remove and discard the cleared solution from each well. Do not touch the beads while removing the solution.



## Step 3. Purify the sample using AMPure XP beads

- 8** Continue to keep the plate or strip tube in the magnetic stand while you dispense 200  $\mu$ l of freshly-prepared 70% ethanol in each sample well.
- 9** Wait for 1 minute to allow any disturbed beads to settle, then remove the ethanol.
- 10** Repeat [step 8](#) to [step 9](#) once.
- 11** Seal the wells with strip caps, then briefly spin the samples to collect the residual ethanol. Return the plate or strip tube to the magnetic stand for 30 seconds. Remove the residual ethanol with a P20 pipette.
- 12** Dry the samples by placing the unsealed plate or strip tube on the thermal cycler, set to hold samples at 37°C, until the residual ethanol has just evaporated (typically 1–2 minutes).

**NOTE**

Do not dry the bead pellet to the point that the pellet appears cracked during any of the bead drying steps in the protocol. Elution efficiency is significantly decreased when the bead pellet is excessively dried.

---

- 13** Add 35  $\mu$ l nuclease-free water to each sample well.
- 14** Seal the wells with strip caps, then mix well on a vortex mixer and briefly spin the plate or strip tube to collect the liquid.
- 15** Incubate for 2 minutes at room temperature.
- 16** Put the plate or strip tube in the magnetic stand and leave for approximately 5 minutes, until the solution is clear.
- 17** Remove the cleared supernatant (approximately 34.5  $\mu$ l) to a fresh PCR plate or strip tube sample well and keep on ice. You can discard the beads at this time.

**NOTE**

It may not be possible to recover the entire 34.5- $\mu$ l supernatant volume at this step; transfer the maximum possible amount of supernatant for further processing. To maximize recovery, transfer the cleared supernatant to a fresh well in two rounds of pipetting, using a P20 pipette set at 17.25  $\mu$ l.

---

### 3 Library Preparation

#### Step 4. Amplify the adaptor-ligated library

## Step 4. Amplify the adaptor-ligated library

This step uses the components listed in [Table 20](#). Before you begin, thaw the reagents listed below and keep on ice. Before use, mix each component as directed.

**Table 20** Reagents for pre-capture PCR amplification

Component	Storage Location	Mixing Method	Where Used
Herculase II Fusion DNA Polymerase (red cap)	SureSelect XT HS and XT Low Input Library Preparation Kit for ILM (Pre PCR), –20°C	Pipette up and down 15–20 times	<a href="#">page 44</a>
5× Herculase II Reaction Buffer (clear cap)	SureSelect XT HS and XT Low Input Library Preparation Kit for ILM (Pre PCR), –20°C	Vortexing	<a href="#">page 44</a>
100 mM dNTP Mix (green cap)	SureSelect XT HS and XT Low Input Library Preparation Kit for ILM (Pre PCR), –20°C	Vortexing	<a href="#">page 44</a>
Forward Primer (brown cap)	SureSelect XT HS and XT Low Input Library Preparation Kit for ILM (Pre PCR), –20°C	Vortexing	<a href="#">page 44</a>
SureSelect XT Low Input Index Primers (yellow plate)	SureSelect XT Low Input Index Primers for ILM (Pre PCR), –20°C	Vortexing	<a href="#">page 44</a>

## Step 4. Amplify the adaptor-ligated library

- 1 Preprogram a thermal cycler (with the heated lid ON) with the program in [Table 21](#). Immediately pause the program, and keep paused until samples are loaded in [step 5](#).

**Table 21** Pre-Capture PCR Thermal Cycler Program\*

Segment	Number of Cycles	Temperature	Time
1	1	98°C	2 minutes
2	8 to 14, based on input DNA quality and quantity (see <a href="#">Table 22</a> )	98°C	30 seconds
		60°C	30 seconds
		72°C	1 minute
3	1	72°C	5 minutes
4	1	4°C	Hold

\* Use a reaction volume setting of 50 µl, if required for thermal cycler set up.

**Table 22** Pre-capture PCR cycle number recommendations

Quality of Input DNA	Quantity of Input DNA	Cycles
Intact DNA from fresh sample	100 to 200 ng	8 cycles
	50 ng	9 cycles
	10 ng	11 cycles
FFPE sample DNA	100 to 200 ng*	11 cycles
	50 ng*	12 cycles
	10 ng*	14 cycles

\* qPCR-determined quantity of amplifiable DNA or DIN value-adjusted amount of input DNA

**CAUTION**

To avoid cross-contaminating libraries, set up PCR reactions (all components except the library DNA) in a dedicated clean area or PCR hood with UV sterilization and positive air flow.

### 3 Library Preparation

#### Step 4. Amplify the adaptor-ligated library

- 2 Prepare the appropriate volume of pre-capture PCR reaction mix, as described in [Table 23](#), on ice. Mix well on a vortex mixer.

**Table 23** Preparation of Pre-Capture PCR Reaction Mix

Reagent	Volume for 1 reaction	Volume for 8 reactions (includes excess)	Volume for 24 reactions (includes excess)
5× Herculase II Reaction Buffer (clear cap)	10 µl	90 µl	250 µl
100 mM dNTP Mix (green cap)	0.5 µl	4.5 µl	12.5 µl
Forward Primer (brown cap)	2 µl	18 µl	50 µl
Herculase II Fusion DNA Polymerase (red cap)	1 µl	9 µl	25 µl
<b>Total</b>	<b>13.5 µl</b>	<b>121.5 µl</b>	<b>337.5 µl</b>

- 3 Add 13.5 µl of the PCR reaction mixture prepared in [Table 23](#) to each purified DNA library sample (34.5 µl) in the PCR plate wells.
- 4 Add 2 µl of the appropriate SureSelect XT Low Input Index Primer (yellow plate; containing P7 index) to each reaction. Use the same P7 index number/plate position as the index number/plate position of the specific P5 Indexed Adaptor ligated to the library in [step 4](#) on [page 39](#). Cap the wells then vortex at high speed for 5 seconds. Spin the plate or strip tube briefly to collect the liquid release any bubbles.
- 5 Before adding the samples to the thermal cycler, resume the program in [Table 21](#) to bring the temperature of the thermal block to 98°C. Once the cycler has reached 98°C, immediately place the sample plate or strip tube in the thermal block and close the lid.

#### CAUTION

The lid of the thermal cycler is hot and can cause burns. Use caution when working near the lid.

## Step 5. Purify the amplified library with AMPure XP beads

In this step, the amplified DNA libraries are purified using AmpPure XP Beads. Critical purification protocol parameters are summarized for experienced users in [Table 24](#).

**Table 24** AMPure XP bead cleanup parameters after pre-capture PCR

Parameter	Value
Volume of RT AMPure XP bead suspension added to each sample well	50 $\mu$ l
Final elution solvent and volume	15 $\mu$ l nuclease-free water
Amount of eluted sample transferred to fresh well	Approximately 15 $\mu$ l

- 1 Verify that the AMPure XP beads were held at room temperature for at least 30 minutes before use.
- 2 Prepare 400  $\mu$ l of 70% ethanol per sample, plus excess, for use in [step 8](#).
- 3 Mix the AMPure XP bead suspension well so that the reagent appears homogeneous and consistent in color.
- 4 Add 50  $\mu$ l of homogeneous AMPure XP beads to each 50- $\mu$ l amplification reaction in the PCR plate or strip tube. Pipette up and down 15–20 times or cap the wells and vortex at high speed for 5–10 seconds to mix.
- 5 Incubate samples for 5 minutes at room temperature.
- 6 Put the plate or strip tube into a magnetic separation device. Wait for the solution to clear (approximately 5 minutes).
- 7 Keep the plate or strip tube in the magnetic stand. Carefully remove and discard the cleared solution from each well. Do not touch the beads while removing the solution.
- 8 Continue to keep the plate or strip tube in the magnetic stand while you dispense 200  $\mu$ l of freshly-prepared 70% ethanol into each sample well.
- 9 Wait for 1 minute to allow any disturbed beads to settle, then remove the ethanol.
- 10 Repeat [step 8](#) and [step 9](#) step once.

### 3 Library Preparation

#### Step 5. Purify the amplified library with AMPure XP beads

- 11** Seal the wells with strip caps, then briefly spin the samples to collect the residual ethanol. Return the plate or strip tube to the magnetic stand for 30 seconds. Remove the residual ethanol with a P20 pipette.
- 12** Dry the samples by placing the unsealed plate or strip tube on the thermal cycler, set to hold samples at 37°C, until the residual ethanol has just evaporated (typically 1–2 minutes).
- 13** Add 15 µl nuclease-free water to each sample well.
- 14** Seal the wells with strip caps, then mix well on a vortex mixer and briefly spin the plate or strip tube to collect the liquid.
- 15** Incubate for 2 minutes at room temperature.
- 16** Put the plate or strip tube in the magnetic stand and leave for 2 to 3 minutes, until the solution is clear.
- 17** Remove the cleared supernatant (approximately 15 µl) to a fresh PCR plate or strip tube sample well and keep on ice. You can discard the beads at this time.

#### NOTE

It may not be possible to recover the entire 15-µl supernatant volume at this step; transfer the maximum possible amount of supernatant for further processing.

---

## Step 6. Assess quality and quantity

Sample analysis can be done with either the 2100 Bioanalyzer instrument or an Agilent TapeStation instrument.

### NOTE

Using either analysis method, observation of a low molecular weight peak, in addition to the expected library fragment peak, indicates the presence of adaptor-dimers in the library. Adaptor-dimer removal is not required for libraries that will be target-enriched in later steps of the workflow. However, for libraries being prepared for whole-genome sequencing (not specifically supported by this user guide), samples with an adaptor-dimer peak must be subjected to an additional round of SPRI-purification. To complete, first dilute the sample to 50 µl with nuclease free water, then follow the SPRI purification procedure on [page 45](#).

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

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

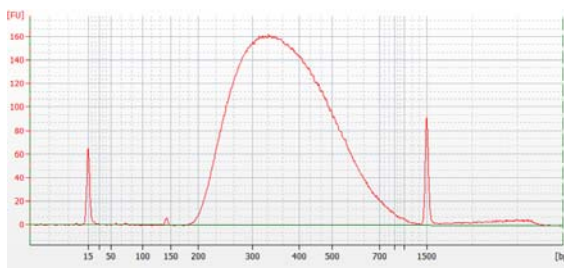
- 1 Set up the 2100 Bioanalyzer instrument as instructed in the reagent kit guide.
- 2 Prepare the chip, samples and ladder as instructed in the reagent kit guide, using 1 µl of each sample for the analysis. Load the prepared chip into the instrument and start the run within five minutes after preparation.
- 3 Verify that the electropherogram shows the peak of DNA fragment size positioned between 300 to 400 bp for high-quality DNA and approximately 200 to 400 bp for FFPE DNA. Sample electropherograms are shown in [Figure 2](#) (library prepared from high-quality DNA), [Figure 3](#) (library prepared from medium-quality FFPE DNA), and [Figure 4](#) (library prepared from low-quality FFPE DNA).

The appearance of an additional low molecular weight peak indicates the presence of adaptor-dimers in the library. It is acceptable to proceed to target enrichment with library samples for which adaptor-dimers are observed in the electropherogram at low abundance, similar to that seen in sample electropherograms on [page 48](#). See Troubleshooting information on [page 103](#) for additional considerations.

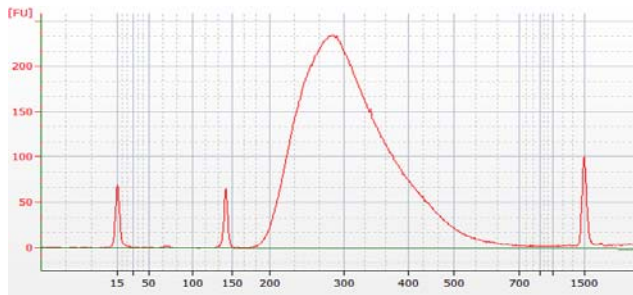
- 4 Determine the concentration of each library by integrating under the entire peak. For accurate quantification, make sure that the concentration falls within the linear range of the assay.

### 3 Library Preparation

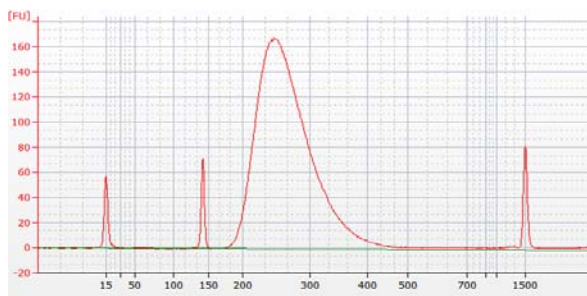
#### Step 6. Assess quality and quantity



**Figure 2** Pre-capture library prepared from a high-quality gDNA sample analyzed using a DNA 1000 Bioanalyzer assay.



**Figure 3** Pre-capture library prepared from a typical FFPE gDNA sample analyzed using a DNA 1000 Bioanalyzer assay.



**Figure 4** Pre-capture library prepared from a low-quality FFPE gDNA sample analyzed using a DNA 1000 Bioanalyzer assay.

**Stopping Point** If you do not continue to the next step, seal the sample wells and store at 4°C overnight or at -20°C for prolonged storage.



**Option 2: Analysis using an Agilent TapeStation and D1000 ScreenTape**

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

- 1 Prepare the TapeStation samples as instructed in the instrument user manual. Use 1  $\mu$ l of each DNA sample diluted with 3  $\mu$ l of D1000 sample buffer for the analysis.

**CAUTION**

For accurate quantitation, make sure to thoroughly mix the combined DNA and sample buffer by vortexing the assay plate or tube strip for 1 minute on the IKA MS3 vortex mixer provided with the 4200/4150 TapeStation system before loading the samples.

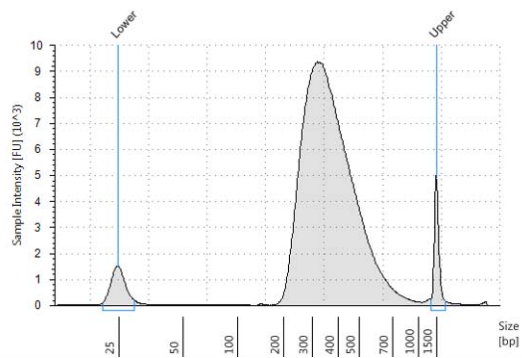
- 2 Load the sample plate or tube strips from [step 1](#), the D1000 ScreenTape, and loading tips into the TapeStation as instructed in the instrument user manual. Start the run.
- 3 Verify that the electropherogram shows the peak of DNA fragment size positioned between 300 to 400 bp for high-quality DNA and approximately 200 to 400 bp for FFPE DNA. Sample electropherograms are shown in [Figure 5](#) (library prepared from high-quality DNA), [Figure 6](#) (library prepared from medium-quality FFPE DNA), and [Figure 7](#) (library prepared from low-quality FFPE DNA).

The appearance of an additional low molecular weight peak indicates the presence of adaptor-dimers in the library. It is acceptable to proceed to target enrichment with library samples for which adaptor-dimers are observed in the electropherogram at low abundance, similar to that seen in sample electropherograms on [page 50](#) to [page 51](#). See Troubleshooting information on [page 103](#) for additional considerations.

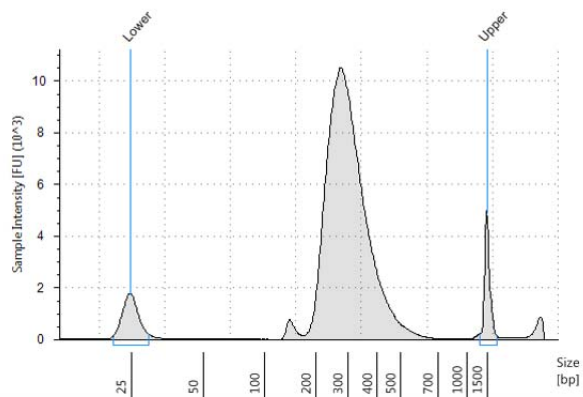
- 4 Determine the concentration of the library DNA by integrating under the peak.

### 3 Library Preparation

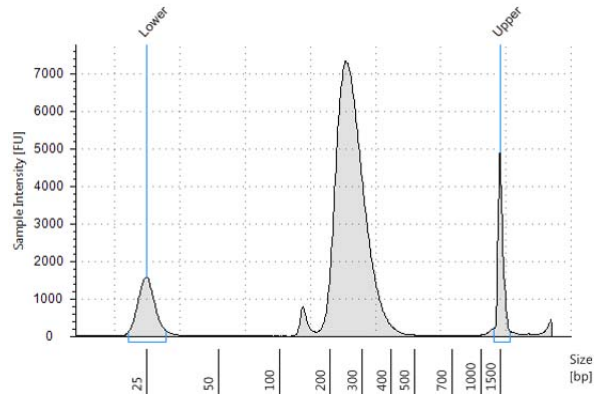
#### Step 6. Assess quality and quantity



**Figure 5** Pre-capture library prepared from a high-quality gDNA sample analyzed using a D1000 ScreenTape assay.



**Figure 6** Pre-capture library prepared from a typical FFPE gDNA sample analyzed using a D1000 ScreenTape assay.



**Figure 7** Pre-capture library prepared from a low-quality FFPE gDNA sample analyzed using a D1000 ScreenTape assay.

**Stopping Point** If you do not continue to the next step, seal the sample wells and store at 4°C overnight or at -20°C for prolonged storage.

### **3 Library Preparation**

Step 6. Assess quality and quantity



## 4 Hybridization and Capture

- Step 1. Hybridize DNA samples to the probe 54
- Step 2. Prepare streptavidin-coated magnetic beads 59
- Step 3. Capture the hybridized DNA using streptavidin-coated beads 60

This chapter describes the steps to hybridize the prepared gDNA libraries with a target-specific probe. After hybridization, the targeted molecules are captured on streptavidin beads. Each DNA library sample is hybridized and captured individually.

The standard single-day protocol includes the hybridization step (approximately 90 minutes) immediately followed by capture and amplification steps. If required, the hybridized samples may be held overnight with capture and amplification steps completed the following day by using the simple protocol modifications noted on [page 55](#).

### CAUTION

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



## 4 Hybridization and Capture

### Step 1. Hybridize DNA samples to the probe

## Step 1. Hybridize DNA samples to the probe

In this step, the prepared gDNA libraries are hybridized to a target-specific probe. For each sample library prepared, do one hybridization and capture. Do not pool samples at this stage.

The hybridization reaction requires 500–1000 ng of prepared DNA in a volume of 12  $\mu$ l. Use the maximum amount of prepared DNA available within this range.

This step uses the components listed in [Table 25](#). Thaw each component under the conditions indicated in the table. Vortex each reagent to mix, then spin tubes briefly to collect the liquid.

**Table 25** Reagents for Hybridization

Kit Component	Storage Location	Thawing Conditions	Where Used
SureSelect XT HS and XT Low Input Blocker Mix (blue cap)	SureSelect XT HS and XT Low Input Target Enrichment Kit ILM Hyb Module, Box 2 (Post PCR), $-20^{\circ}\text{C}$	Thaw on ice	<a href="#">page 55</a>
SureSelect RNase Block (purple cap)	SureSelect XT HS and XT Low Input Target Enrichment Kit ILM Hyb Module, Box 2 (Post PCR), $-20^{\circ}\text{C}$	Thaw on ice	<a href="#">page 56</a>
SureSelect Fast Hybridization Buffer (bottle)	SureSelect XT HS and XT Low Input Target Enrichment Kit ILM Hyb Module, Box 2 (Post PCR), $-20^{\circ}\text{C}$	Thaw and keep at Room Temperature	<a href="#">page 57</a>
Probe	$-80^{\circ}\text{C}$	Thaw on ice	<a href="#">page 57</a>

- 1 Preprogram a thermal cycler (with the heated lid ON) with the program in [Table 26](#). Immediately pause the program, and keep paused until samples are loaded in [step 4](#) on [page 56](#).

**Table 26** Pre-programmed thermal cycler program for Hybridization\*

Segment #	Number of Cycles	Temperature	Time
1	1	95°C	5 minutes
2	1	65°C	10 minutes
3	1	65°C	1 minute (Pause cycler here for reagent addition; see <a href="#">step 7</a> on <a href="#">page 58</a> )
4	60	65°C <sup>†</sup>	1 minute
		37°C	3 seconds
5	1	65°C <sup>†</sup>	Hold briefly until ready to begin capture steps on <a href="#">page 60</a>

\* Use a reaction volume setting of 30 µl (final volume of hybridization reactions during cycling in Segment 4).

† Hybridization at 65°C is optimal for probes designed for the SureSelect XT HS2/XT HS/XT Low Input platforms. Reducing the hybridization temperature (Segments 4 and 5) may improve performance for probes designed for the SureSelect XT platform, including SureSelect XT Human All Exon V6 (62.5°C), SureSelect XT Clinical Research Exome V2 (62.5°C) and custom probes originally designed for use with SureSelect XT system (60°C–65°C).

**NOTE**

The Hybridization thermal cycling program in [Table 26](#) requires about 90 minutes. The Hybridization reaction may be run overnight with the following protocol modifications:

- In segment 5 of the thermal cycler program ([Table 26](#)), replace the 65°C Hold step with a 21°C Hold step.
- The hybridized samples may be held at 21°C for up to 16 hours. Begin the capture preparation steps on [page 59](#) on day 2, after the overnight hold.

- 2 Place 500–1000 ng of each prepared gDNA library sample into the hybridization plate or strip tube wells and then bring the final volume in each well to 12 µl using nuclease-free water. Use the maximum possible amount of each prepped DNA, within the 500–1000 ng range.
- 3 To each DNA library sample well, add 5 µl of SureSelect XT HS and XT Low Input Blocker Mix (blue cap). Seal the wells then vortex at high speed for 5 seconds. Spin the plate or strip tube briefly to collect the liquid and release any bubbles.

## 4 Hybridization and Capture

### Step 1. Hybridize DNA samples to the probe

#### CAUTION

The lid of the thermal cycler is hot and can cause burns. Use caution when working near the lid.

- 4 Transfer the sealed sample plate or strip to the thermal cycler and resume the thermal cycling program (Table 26 on page 55), allowing the cycler to complete Segments 1 and 2 of the program.

**Important: The thermal cycler must be paused during Segment 3 to allow additional reagents to be added to the Hybridization wells in step 7 on page 58.**

During Segments 1 and 2 of the thermal cycling program, begin preparing the additional hybridization reagents as described in step 5 and step 6 below. If needed, you can finish these preparation steps after pausing the thermal cycler in Segment 3.

- 5 Prepare a 25% solution of SureSelect RNase Block (1 part RNase Block to 3 parts water) according to Table 27. Prepare the amount required for the number of hybridization reactions in the run, plus excess. Mix well and keep on ice.

**Table 27** Preparation of RNase Block solution

Reagent	Volume for 1 reaction	Volume for 8 reactions (includes excess)	Volume for 24 reactions (includes excess)
SureSelect RNase Block (purple cap)	0.5 µl	4.5 µl	12.5 µl
Nuclease-free water	1.5 µl	13.5 µl	37.5 µl
<b>Total</b>	<b>2 µl</b>	<b>18 µl</b>	<b>50 µl</b>



**NOTE**

Prepare the mixture described in [step 6](#), below, just before pausing the thermal cycler in Segment 3. Keep the mixture at room temperature briefly until the mixture is added to the DNA samples in [step 7](#) on [page 58](#). Do not keep solutions containing the probe at room temperature for extended periods.

- 6** Prepare the Probe Hybridization Mix appropriate for your probe design size. Use [Table 28](#) for probes  $\geq 3$  Mb or [Table 29](#) for probes  $< 3$  Mb.

**Combine the listed reagents at room temperature.** Mix well by vortexing at high speed for 5 seconds then spin down briefly. Proceed immediately to [step 7](#).

**Table 28** Preparation of Probe Hybridization Mix for **probes  $\geq 3$  Mb**

Reagent	Volume for 1 reaction	Volume for 8 reactions (includes excess)	Volume for 24 reactions (includes excess)
25% RNase Block solution (from <a href="#">step 5</a> )	2 $\mu$ l	18 $\mu$ l	50 $\mu$ l
Probe (with design $\geq 3$ Mb)	5 $\mu$ l	45 $\mu$ l	125 $\mu$ l
SureSelect Fast Hybridization Buffer	6 $\mu$ l	54 $\mu$ l	150 $\mu$ l
<b>Total</b>	<b>13 <math>\mu</math>l</b>	<b>117 <math>\mu</math>l</b>	<b>325 <math>\mu</math>l</b>

**Table 29** Preparation of Probe Hybridization Mix for **probes  $< 3$  Mb**

Reagent	Volume for 1 reaction	Volume for 8 reactions (includes excess)	Volume for 24 reactions (includes excess)
25% RNase Block solution (from <a href="#">step 5</a> )	2 $\mu$ l	18 $\mu$ l	50 $\mu$ l
Probe (with design $< 3$ Mb)	2 $\mu$ l	18 $\mu$ l	50 $\mu$ l
SureSelect Fast Hybridization Buffer	6 $\mu$ l	54 $\mu$ l	150 $\mu$ l
Nuclease-free water	3 $\mu$ l	27 $\mu$ l	75 $\mu$ l
<b>Total</b>	<b>13 <math>\mu</math>l</b>	<b>117 <math>\mu</math>l</b>	<b>325 <math>\mu</math>l</b>

## 4 Hybridization and Capture

### Step 1. Hybridize DNA samples to the probe

- 7 Once the thermal cycler starts Segment 3 of the program in [Table 26](#) (1 minute at 65°C), pause the program. With the cycler paused, and while keeping the DNA + Blocker samples in the cycler, transfer 13 µl of the room-temperature Probe Hybridization Mix from [step 6](#) to each sample well.

Mix well by pipetting up and down slowly 8 to 10 times.

The hybridization reaction wells now contain approximately 30 µl.

- 8 Seal the wells with fresh strip caps. Make sure that all wells are completely sealed. Vortex briefly, then spin the plate or strip tube briefly to remove any bubbles from the bottom of the wells. Immediately return the plate or strip tube to the thermal cycler.
- 9 Resume the thermal cycling program to allow hybridization of the prepared DNA samples to the Probe.

#### CAUTION

Wells must be adequately sealed to minimize evaporation, or your results can be negatively impacted.

Before you do the first experiment, make sure the plasticware and capping method are appropriate for the thermal cycler. Check that no more than 4 µl is lost to evaporation under the conditions used for hybridization.

---

## Step 2. Prepare streptavidin-coated magnetic beads

The remaining hybridization capture steps use the reagents in [Table 30](#).

### NOTE

If performing same-day hybridization and capture, begin the bead preparation steps below approximately one hour after starting hybridization in [step 9](#) on [page 58](#). If performing next-day capture after an overnight hold at 21°C, begin the bead preparation steps below on day 2, just before you are ready to start the capture steps on [page 60](#).

**Table 30** Reagents for Capture

Kit Component	Storage Location	Where Used
SureSelect Binding Buffer	SureSelect Target Enrichment Kit, ILM Hyb Module, Box 1 (Post PCR), RT	<a href="#">page 59</a>
SureSelect Wash Buffer 1	SureSelect Target Enrichment Kit, ILM Hyb Module, Box 1 (Post PCR), RT	<a href="#">page 60</a>
SureSelect Wash Buffer 2	SureSelect Target Enrichment Kit, ILM Hyb Module, Box 1 (Post PCR), RT	<a href="#">page 60</a>
Dynabeads MyOne Streptavidin T1	Follow storage recommendations provided by supplier (see <a href="#">Table 4</a> on <a href="#">page 15</a> )	<a href="#">page 59</a>

- 1 Vigorously resuspend the Dynabeads MyOne Streptavidin T1 magnetic beads on a vortex mixer. The magnetic beads settle during storage.
- 2 For each hybridization sample, add 50 µl of the resuspended beads to wells of a fresh PCR plate or a strip tube.
- 3 Wash the beads:
  - a Add 200 µl of SureSelect Binding Buffer.
  - b Mix by pipetting up and down 20 times or cap the wells and vortex at high speed for 5–10 seconds then spin down briefly.
  - c Put the plate or strip tube into a magnetic separator device.
  - d Wait at least 5 minutes or until the solution is clear, then remove and discard the supernatant.
  - e Repeat [step a](#) through [step d](#) two more times for a total of 3 washes.
- 4 Resuspend the beads in 200 µl of SureSelect Binding Buffer.

### NOTE

If you are equipped for higher-volume magnetic bead captures, the streptavidin beads may instead be batch-washed in an Eppendorf tube or conical vial.

## 4 Hybridization and Capture

### Step 3. Capture the hybridized DNA using streptavidin-coated beads

## Step 3. Capture the hybridized DNA using streptavidin-coated beads

- 1 After all streptavidin bead preparation steps are complete, and with the hybridization thermal cycling program in the final hold segment (see [Table 26](#) on page 55), transfer the samples to room temperature.
- 2 Immediately transfer the entire volume (approximately 30  $\mu$ l) of each hybridization mixture to wells containing 200  $\mu$ l of washed streptavidin beads using a multichannel pipette.  
Pipette up and down 5–8 times to mix then seal the wells with fresh caps.
- 3 Incubate the capture plate or strip tube on a 96-well plate mixer, mixing vigorously (at 1400–1800 rpm), for 30 minutes at room temperature.  
**Make sure the samples are properly mixing in the wells.**
- 4 During the 30-minute incubation for capture, prewarm SureSelect Wash Buffer 2 at 70°C as described below.
  - a Place 200- $\mu$ l aliquots of Wash Buffer 2 in wells of a fresh 96-well plate or strip tubes. Aliquot 6 wells of buffer for each DNA sample in the run.
  - b Cap the wells and then incubate in the thermal cycler, with heated lid ON, held at 70°C until used in [step 9](#).
- 5 When the 30-minute incubation period initiated in [step 3](#) is complete, spin the samples briefly to collect the liquid.
- 6 Put the plate or strip tube in a magnetic separator to collect the beads. Wait until the solution is clear, then remove and discard all of the supernatant.
- 7 Resuspend the beads in 200  $\mu$ l of SureSelect Wash Buffer 1. Mix by pipetting up and down 15–20 times, until beads are fully resuspended.
- 8 Put the plate or strip tube in the magnetic separator. Wait for the solution to clear (approximately 1 minute), then remove and discard all of the supernatant.

## Step 3. Capture the hybridized DNA using streptavidin-coated beads

**CAUTION**

It is important to maintain bead suspensions at 70°C during the washing procedure below to ensure specificity of capture.

Make sure that the SureSelect Wash Buffer 2 is pre-warmed to 70°C before use.

Do not use a tissue incubator, or other devices with significant temperature fluctuations, for the incubation steps.

- 
- 9 Remove the plate or strip tubes from the magnetic separator and transfer to a rack at room temperature.** Wash the beads with Wash Buffer 2, using the protocol steps below.
- a** Resuspend the beads in 200 µl of 70°C prewarmed Wash Buffer 2. Pipette up and down 15–20 times, until beads are fully resuspended.
  - b** Seal the wells with fresh caps and then vortex at high speed for 8 seconds. Spin the plate or strip tube briefly to collect the liquid without pelleting the beads.  
**Make sure the beads are in suspension before proceeding.**
  - c** Incubate the samples for 5 minutes at 70°C in the thermal cycler with the heated lid ON.
  - d** Put the plate or strip tube in the magnetic separator at room temperature.
  - e** Wait 1 minute for the solution to clear, then remove and discard the supernatant.
  - f** Repeat [step a](#) through [step e](#) five more times for a total of 6 washes.
- 10** After verifying that all wash buffer has been removed, add 25 µl of nuclease-free water to each sample well. Pipette up and down 8 times to resuspend the beads.

Keep the samples on ice until they are used on [page 66](#).

**NOTE**

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

## **4 Hybridization and Capture**

**Step 3. Capture the hybridized DNA using streptavidin-coated beads**



## 5 Post-Capture Sample Processing for Multiplexed Sequencing

- Step 1. Amplify the captured libraries 64
- Step 2. Purify the amplified captured libraries using AMPure XP beads 67
- Step 3. Assess sequencing library DNA quantity and quality 69
- Step 4. Pool samples for multiplexed sequencing 74
- Step 5. Prepare sequencing samples 76
- Step 6. Do the sequencing run and analyze the data 78
- Sequence analysis resources 83

This chapter describes the steps to amplify, purify, and assess quality and quantity of the captured libraries. Sample pooling instructions are provided to prepare the dual-indexed samples for multiplexed sequencing.



## 5 Post-Capture Sample Processing for Multiplexed Sequencing

### Step 1. Amplify the captured libraries

## Step 1. Amplify the captured libraries

In this step, the SureSelect-enriched DNA libraries are PCR amplified.

This step uses the components listed in [Table 31](#). Before you begin, thaw the reagents listed below and keep on ice.

**Table 31** Reagents for post-capture PCR amplification

Component	Storage Location	Mixing Method	Where Used
Herculase II Fusion DNA Polymerase (red cap)	SureSelect XT HS and XT Low Input Target Enrichment Kit, ILM Hyb Module Box 2 (Post PCR), -20°C	Pipette up and down 15–20 times	<a href="#">page 66</a>
5× Herculase II Reaction Buffer (clear cap)	SureSelect XT HS and XT Low Input Target Enrichment Kit, ILM Hyb Module Box 2 (Post PCR), -20°C	Vortexing	<a href="#">page 66</a>
100 mM dNTP Mix (green cap)	SureSelect XT HS and XT Low Input Target Enrichment Kit, ILM Hyb Module Box 2 (Post PCR), -20°C	Vortexing	<a href="#">page 66</a>
SureSelect Post-Capture Primer Mix (clear cap)	SureSelect XT HS and XT Low Input Target Enrichment Kit, ILM Hyb Module Box 2 (Post PCR), -20°C	Vortexing	<a href="#">page 66</a>

Prepare one amplification reaction for each DNA library.

### CAUTION

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



- 1 Preprogram a thermal cycler (with the heated lid ON) with the program in [Table 32](#). Immediately pause the program, and keep paused until samples are loaded in [step 5](#).

**Table 32** Post-capture PCR Thermal Cycler Program

Segment	Number of Cycles	Temperature	Time
1	1	98°C	2 minutes
2	9 to 14 See <a href="#">Table 33</a> for hybridization probe design size-based cycle number recommendations	98°C	30 seconds
		60°C	30 seconds
		72°C	1 minute
3	1	72°C	5 minutes
4	1	4°C	Hold

**Table 33** Post-capture PCR cycle number recommendations

Probe Size/Description	Cycles
Probes <0.2 Mb	14 cycles
Probes 0.2–3 Mb (includes ClearSeq Comp Cancer)	12 cycles
Probes 3–5 Mb	10 cycles
Probes >5 Mb (includes Human All Exon probes)	9 cycles

## 5 Post-Capture Sample Processing for Multiplexed Sequencing

### Step 1. Amplify the captured libraries

- 2 Prepare the appropriate volume of PCR reaction mix, as described in [Table 34](#), on ice. Mix well on a vortex mixer.

**Table 34** Preparation of post-capture PCR Reaction mix

Reagent	Volume for 1 reaction	Volume for 8 reactions (includes excess)	Volume for 24 reactions (includes excess)
Nuclease-free water	12.5 µl	112.5 µl	312.5 µl
5× Herculase II Reaction Buffer (clear cap)	10 µl	90 µl	250 µl
Herculase II Fusion DNA Polymerase (red cap)	1 µl	9 µl	25 µl
100 mM dNTP Mix (green cap)	0.5 µl	4.5 µl	12.5 µl
SureSelect Post-Capture Primer Mix (clear cap)	1 µl	9 µl	25 µl
<b>Total</b>	<b>25 µl</b>	<b>225 µl</b>	<b>625 µl</b>

- 3 Add 25 µl of the PCR reaction mix prepared in [Table 34](#) to each sample well containing 25 µl of bead-bound target-enriched DNA (prepared on [page 61](#) and held on ice).
- 4 Mix the PCR reactions well by pipetting up and down until the bead suspension is homogeneous. Avoid splashing samples onto well walls; do not spin the samples at this step.
- 5 Place the plate or strip tube in the thermal cycler and resume the thermal cycling program in [Table 32](#).
- 6 When the PCR amplification program is complete, spin the plate or strip tube briefly. Remove the streptavidin-coated beads by placing the plate or strip tube on the magnetic stand at room temperature. Wait 2 minutes for the solution to clear, then **remove each supernatant (approximately 50 µl) to wells of a fresh plate or strip tube.**  
The beads can be discarded at this time.

## Step 2. Purify the amplified captured libraries using AMPure XP beads

In this step, the amplified enriched DNA libraries are purified using AmpPure XP Beads. Critical purification protocol parameters are summarized for experienced users in [Table 35](#).

**Table 35** AMPure XP bead cleanup parameters after post-capture PCR

Parameter	Value
Volume of RT AMPure XP bead suspension added to each sample well	50 $\mu$ l
Final elution solvent and volume	25 $\mu$ l nuclease-free water
Amount of eluted sample transferred to fresh well	Approximately 25 $\mu$ l

- 1 Let the AMPure XP beads come to room temperature for at least 30 minutes.
- 2 Prepare 400  $\mu$ l of fresh 70% ethanol per sample, plus excess, for use in [step 8](#).
- 3 Mix the AMPure XP bead suspension well so that the suspension appears homogeneous and consistent in color.
- 4 Add 50  $\mu$ l of the homogeneous AMPure XP bead suspension to each amplified DNA sample (approximately 50  $\mu$ l) in the PCR plate or strip tube. Mix well by pipetting up and down 15–20 times or cap the wells and vortex at high speed for 5–10 seconds.  
 Check that the beads are in a homogeneous suspension in the sample wells. Each well should have a uniform color with no layers of beads or clear liquid present.
- 5 Incubate samples for 5 minutes at room temperature.
- 6 Put the plate or strip tube on the magnetic stand at room temperature. Wait for the solution to clear (approximately 3 to 5 minutes).
- 7 While keeping the plate or tubes in the magnetic stand, carefully remove and discard the cleared solution from each well. Do not disturb the beads while removing the solution.
- 8 Continue to keep the plate or tubes in the magnetic stand while you dispense 200  $\mu$ l of freshly-prepared 70% ethanol in each sample well.

## 5 Post-Capture Sample Processing for Multiplexed Sequencing

### Step 2. Purify the amplified captured libraries using AMPure XP beads

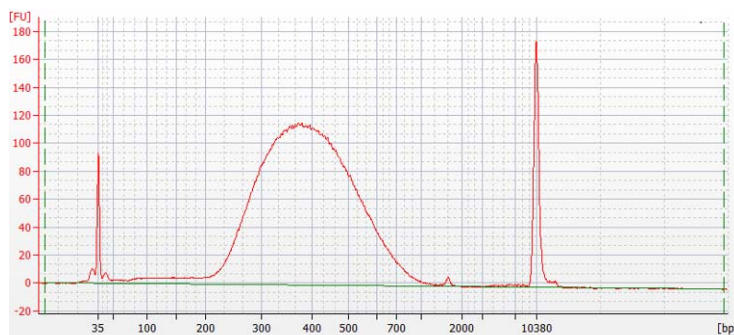
- 9 Wait for 1 minute to allow any disturbed beads to settle, then remove the ethanol.
- 10 Repeat [step 8](#) and [step 9](#) once for a total of two washes. Make sure to remove all of the ethanol at each wash step.
- 11 Seal the wells with strip caps, then briefly spin to collect the residual ethanol. Return the plate or strip tube to the magnetic stand for 30 seconds. Remove the residual ethanol with a P20 pipette.
- 12 Dry the samples by placing the unsealed plate or strip tube on the thermal cycler, set to hold samples at 37°C, until the residual ethanol has just evaporated (typically 1–2 minutes).
- 13 Add 25 µl of nuclease-free water to each sample well.
- 14 Seal the sample wells, then mix well on a vortex mixer and briefly spin to collect the liquid without pelleting the beads.
- 15 Incubate for 2 minutes at room temperature.
- 16 Put the plate or strip tube in the magnetic stand and leave for 2 minutes or until the solution is clear.
- 17 Remove the cleared supernatant (approximately 25 µl) to a fresh well. You can discard the beads at this time.

## Step 3. Assess sequencing library DNA quantity and quality

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

Use the Bioanalyzer High Sensitivity DNA Assay to analyze the amplified captured DNA. Perform the assay according to the [High Sensitivity DNA Kit Guide](#).

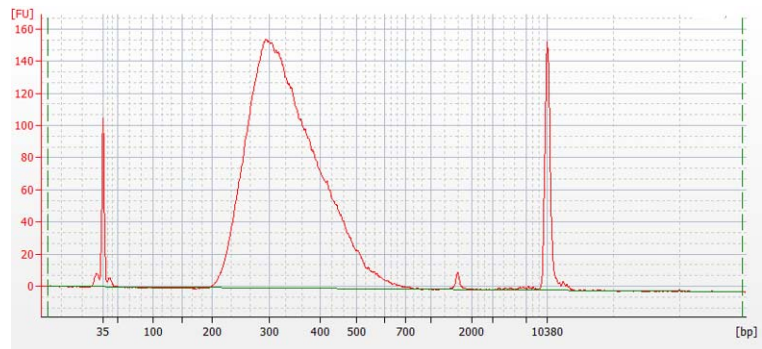
- 1 Set up the 2100 Bioanalyzer instrument as instructed in the reagent kit guide.
- 2 Prepare the chip, samples and ladder as instructed in the reagent kit guide, using 1  $\mu$ l of each sample for the analysis.
- 3 Load the prepared chip into the instrument and start the run within five minutes after preparation.
- 4 Verify that the electropherogram shows the peak of DNA fragment size positioned between 200 and 400 bp. Sample electropherograms are shown in [Figure 8](#) (library prepared from high-quality DNA), [Figure 9](#) (library prepared from medium-quality FFPE DNA), and [Figure 10](#) (library prepared from low-quality FFPE DNA).
- 5 Measure the concentration of each library by integrating under the entire peak. For accurate quantification, make sure that the concentration falls within the linear range of the assay.



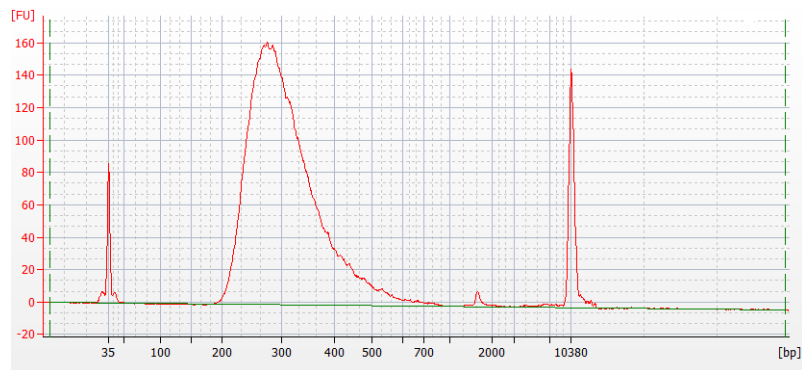
**Figure 8** Post-capture library prepared from a high-quality gDNA sample analyzed using a Bioanalyzer system High Sensitivity DNA assay.

## 5 Post-Capture Sample Processing for Multiplexed Sequencing

### Step 3. Assess sequencing library DNA quantity and quality



**Figure 9** Post-capture library prepared from a typical FFPE gDNA sample analyzed using a Bioanalyzer system High Sensitivity DNA assay.



**Figure 10** Post-capture library prepared from a low-quality FFPE gDNA sample analyzed using a Bioanalyzer system High Sensitivity DNA assay.

**Stopping Point** If you do not continue to the next step, seal the plate and store at 4°C overnight or at -20°C for prolonged storage.

### Option 2: Analysis using an Agilent TapeStation and High Sensitivity D1000 ScreenTape

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

- 1 Prepare the TapeStation samples as instructed in the in the reagent kit guide. Use 2  $\mu$ l of each dual-indexed DNA sample diluted with 2  $\mu$ l of High Sensitivity D1000 sample buffer for the analysis.

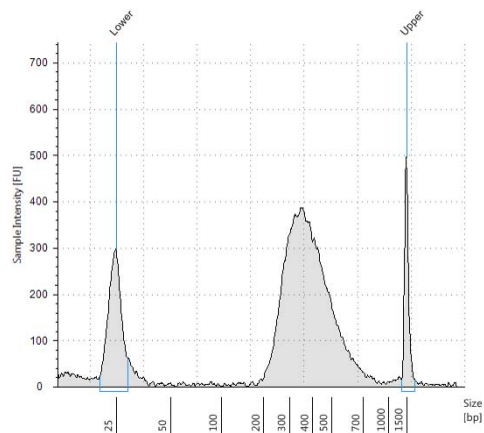
#### CAUTION

For accurate quantitation, make sure to thoroughly mix the combined DNA and sample buffer by vortexing the assay plate or tube strip for 1 minute on the IKA MS3 vortex mixer provided with the 4200/4150 TapeStation system before loading the samples.

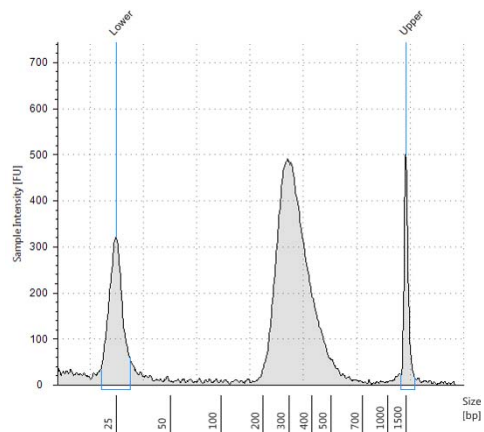
- 2 Load the sample plate or tube strips from [step 1](#), the High Sensitivity D1000 ScreenTape, and loading tips into the TapeStation as instructed in the reagent kit guide. Start the run.
- 3 Verify that the electropherogram shows the peak of DNA fragment size positioned between 200 and 400 bp. Sample electropherograms are shown in [Figure 8](#) (library prepared from high-quality DNA), [Figure 9](#) (library prepared from medium-quality FFPE DNA), and [Figure 10](#) (library prepared from low-quality FFPE DNA).
- 4 Determine the concentration of each library by integrating under the entire peak.

## 5 Post-Capture Sample Processing for Multiplexed Sequencing

### Step 3. Assess sequencing library DNA quantity and quality

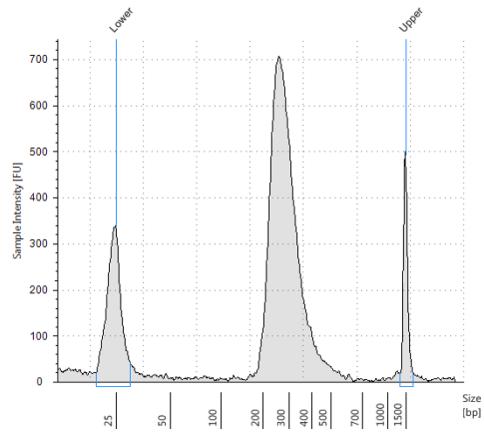


**Figure 11** Post-capture library prepared from a high-quality gDNA sample analyzed using a High Sensitivity D1000 ScreenTape assay.



**Figure 12** Post-capture library prepared from a typical FFPE gDNA sample analyzed using a High Sensitivity D1000 ScreenTape assay.





**Figure 13** Post-capture library prepared from a low-quality FFPE gDNA sample analyzed using a High Sensitivity D1000 ScreenTape assay.

**Stopping Point** If you do not continue to the next step, seal the plate and store at 4°C overnight or at -20°C for prolonged storage.

## Step 4. Pool samples for multiplexed sequencing

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

Combine the libraries such that each dual-indexed sample is present in equimolar amounts in the pool using one of the following methods:

**Method 1:** Dilute each sample to be pooled to the same final concentration (typically 4 nM–15 nM, or the concentration of the most dilute sample) using Low TE, then combine equal volumes of all samples to create the final pool.

**Method 2:** Starting with samples at different concentrations, add the appropriate volume of each sample to achieve equimolar concentration in the pool, then adjust the pool to the desired final volume using Low TE. The formula below is provided for determination of the amount of each sample to add to the pool.

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

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

$C(f)$  is the desired final concentration of all the DNA in the pool (typically 4 nM–15 nM or the concentration of the most dilute sample)

$\#$  is the number of dual-indexed samples, and

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

Table 36 shows an example of the amount of 4 dual-indexed samples (of different concentrations) and Low TE needed for a final volume of 20  $\mu$ l at 10 nM DNA.

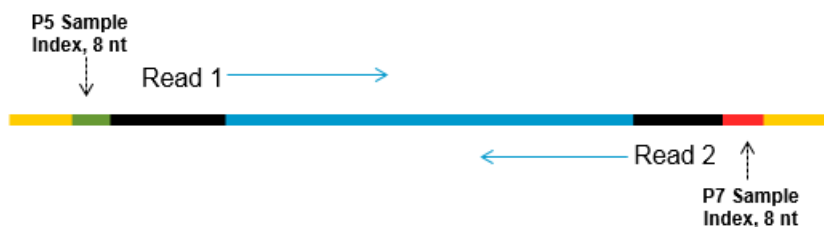
**Table 36** Example of volume calculation for total volume of 20 µl at 10 nM concentration

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

If you store the library before sequencing, add Tween 20 to 0.1% v/v and store at -20°C short term.

## Step 5. Prepare sequencing samples

The final SureSelect<sup>XT</sup> Low Input library pool is ready for direct sequencing using standard Illumina paired-end primers and chemistry. Each fragment in the prepared library contains one target insert surrounded by sequence motifs required for multiplexed sequencing using the Illumina platform, as shown in [Figure 14](#).



**Figure 14** Content of SureSelect XT Low Input dual-indexed sequencing library. Each fragment contains one target insert (blue) surrounded by the Illumina paired-end sequencing elements (black), the P5 and P7 sample indexes (green and red), and the library bridge PCR primers (yellow).

Libraries can be sequenced on the Illumina HiSeq, MiSeq, NextSeq, or NovaSeq platform using the run type and chemistry combinations shown in [Table 37](#).

### CAUTION

Reduced P5 sample index quality has been observed when SureSelect<sup>XT</sup> Low Input libraries are sequenced on the HiSeq2500 instrument in high-output run mode (v4 chemistry). Lower Q scores have been shown to impact coverage and sensitivity of variant calls, especially for aberrations present at less than 10% frequency.

Proceed to cluster amplification using the appropriate Illumina Paired-End Cluster Generation Kit. See [Table 37](#) for kit configurations compatible with the recommended read length.

The optimal seeding concentration for SureSelect<sup>XT</sup> Low Input target-enriched libraries varies according to sequencing platform, run type, and Illumina kit version. See [Table 37](#) for guidelines. Seeding concentration and cluster density may also need to be optimized based on the DNA fragment size range for the library and on the desired output and data quality. Begin optimization using a seeding concentration in the middle of the range listed in [Table 37](#).

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

**Table 37** Illumina Kit Configuration Selection Guidelines

Platform	Run Type	Read Length	SBS Kit Configuration	Chemistry	Seeding Concentration
HiSeq 2500	Rapid Run	2 × 100 bp	200 Cycle Kit	v2	9–10 pM
HiSeq 2500	High Output*	2 × 100 bp	250 Cycle Kit	v4	12–14 pM
MiSeq	All Runs	2 × 100 bp or 2 × 150 bp	300 Cycle Kit	v2	9–10 pM
MiSeq	All Runs	2 × 75 bp	150 Cycle Kit	v3	12–16 pM
NextSeq 500/550	All Runs	2 × 100 bp or 2 × 150 bp	300 Cycle Kit	v2.5	1.2–1.5 pM
HiSeq 3000/4000	All Runs	2 × 100 bp or 2 × 150 bp	300 Cycle Kit	v1	300–400 pM
NovaSeq 6000	Standard Workflow Runs	2 × 100 bp or 2 × 150 bp	300 Cycle Kit	v1.0 or v1.5	300–600 pM
NovaSeq 6000	Xp Workflow Runs	2 × 100 bp or 2 × 150 bp	300 Cycle Kit	v1.0 or v1.5	200–400 pM

\* Reduced i5 index sequence quality and lowered Q scores have been observed in sequences obtained from HiSeq 2500 High Output (v4 chemistry) runs.

#### NOTE

For All-In-One assays that include translocation detection, Agilent strongly recommends using paired-end sequencing read length of at least 2 × 100 bp and preferably 2 × 150 bp.

## Step 6. Do the sequencing run and analyze the data

Use the guidelines below for SureSelect<sup>XT</sup> Low Input dual-indexed library sequencing run setup and analysis.

- Each sample-level index (i7 and i5) requires an 8-bp index read. For complete i7 index sequence information, see [Table 52](#) on page 98. For complete i5 index sequence information, see [Table 53](#) on page 99 and [Table 54](#) on page 100.
- For the HiSeq, NextSeq, and NovaSeq platforms, set up the run using the instrument's user interface, following the guidelines on [page 79](#).
- For the MiSeq platform, set up the run using Illumina Experiment Manager (IEM) using the steps detailed on [page 79](#) to [page 82](#) to generate a custom sample sheet.
- Do not use Illumina's IEM adaptor trimming options. Make sure any IEM adaptor trimming option checkboxes are cleared (deselected) when setting up the sequencing run. Adaptors are trimmed in later processing steps using the Agilent software tools described below to ensure proper processing.
- Demultiplex using Illumina's bcl2fastq, BCL Convert or DRAGEN software to generate paired end reads based on the dual indexes and remove sequences with incorrectly paired P5 and P7 indexes.
- For human germline DNA variant analysis, you can use Agilent's Alissa Reporter software for the complete FASTQ file to variant discovery process (see [page 83](#) for more information).
- For germline or somatic variant analysis, you can use Agilent's AGeNT software modules to process the library read FASTQ files to analysis-ready BAM files. See [page 84](#) for more information.

## HiSeq/NextSeq/NovaSeq platform sequencing run setup guidelines

Set up sequencing runs using the instrument control software interface, using the settings shown in [Table 38](#). For HiSeq runs, select *Dual Index* on the *Run Configuration* screen of the instrument control software interface and enter the **Cycles** settings in [Table 38](#).

For the NextSeq or NovaSeq platform, open the *Run Setup* screen of the instrument control software interface and enter the **Read Length** settings in [Table 38](#). In the **Custom Primers** section, clear (do **not** select) the checkboxes for all primers (*Read 1*, *Read 2*, *Index 1* and *Index 2*).

**Table 38** Run settings

Run Segment	Cycles/Read Length
Read 1	100
Index 1 (i7)	8
Index 2 (i5)	8
Read 2	100

## MiSeq platform sequencing run setup guidelines

Use the Illumina Experiment Manager (IEM) software to generate a custom Sample Sheet according to the guidelines below. Once a Sample Sheet has been generated, index sequences need to be manually changed to the dual indexes used for each sample. See [Table 52](#) on page 98 and [Table 53](#) on page 99 for nucleotide sequences of the dual indexes.

### 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*.

## 5 Post-Capture Sample Processing for Multiplexed Sequencing

### Step 6. Do the sequencing run and analyze the data

- 2 On the **Workflow Parameters** screen, enter the run information, making sure to specify the key parameters highlighted below. In the *Library Prep Workflow* field, select **TruSeq Nano DNA**. In the *Index Adapters* field, select **TruSeq DNA CD Indexes (96 Indexes)**. If your pipeline uses SureCall for adaptor trimming, then make sure to clear both adaptor-trimming checkboxes under *FASTQ Only Workflow-Specific Settings* (circled below), since these are selected by default.

If **TruSeq Nano DNA** is not available in the *Sample Prep Kit* field, instead select **TruSeq HT**.

## Sample Sheet Wizard - Workflow Parameters

FASTQ Only Run Settings

Reagent Cartridge Barcode\* MS5871368-300V2

Library Prep Workflow TruSeq Nano DNA

Index Adapters TruSeq DNA CD Indexes (96 Indexes)

Index Reads  0 (None)  1 (Single)  2 (Dual)

Experiment Name

Investigator Name

Description

Date 1/22/2018

Read Type  Paired End  Single Read

Cycles Read 1 100

Cycles Read 2 100

\* - required field

FASTQ Only Workflow-Specific Settings

Custom Primer for Read 1

Custom Primer for Index

Custom Primer for Read 2

Reverse Complement

Use Adapter Trimming

Use Adapter Trimming Read 2



- 3 Using the **Sample Sheet Wizard**, set up a New Plate, entering the required information for each sample to be sequenced. In the **I7 Sequence** column, assign each sample to any of the Illumina i7 indexes. The index will be corrected to a SureSelect XT Low Input index at a later stage.

Likewise, in the **I5 Sequence** column, assign any of the Illumina i5 indexes, to be corrected to the i5 sequence from the SureSelect XT Low Input Dual Index P5 Indexed Adaptor at a later stage.



EM Illumina Experiment Manager

Illumina Experiment Manager

Sample Sheet Wizard - Sample Selection

Samples to include in sample sheet

Sample ID*	Sample Name	Plate	Well	Index1 (I7)*	I7 Sequence	Index2 (I5)*	I5 Sequence	Sample Project	Description
1	1	Plate1	A01	D701	ATTACTCG	D501	TATAGCCT		
2	2	Plate1	A02	D702	TCCGGAGA	D501	TATAGCCT		
3	3	Plate1	A03	D703	CGCTCATT	D501	TATAGCCT		
4	4	Plate1	A04	D704	GAGATTCC	D501	TATAGCCT		
5	5	Plate1	A05	D705	ATTCAGAA	D501	TATAGCCT		
6	6	Plate1	A06	D706	GAATTCGT	D501	TATAGCCT		

- 4 Finish the sample sheet setup tasks and save the sample sheet file.

## 5 Post-Capture Sample Processing for Multiplexed Sequencing

### Step 6. Do the sequencing run and analyze the data

#### Editing the Sample Sheet to include SureSelect XT Low Input dual indexes

- 1 Open the Sample Sheet file in a text editor and edit the i7 and i5 index information for each sample in columns 5–8 (highlighted below).
  - In column 5 under **I7\_Index\_ID**, enter the name of the SureSelect XT Low Input index assigned to the sample. In column 6 under **index**, enter the corresponding P7 index sequence. See [Table 52](#) on page 98 for nucleotide sequences of the SureSelect XT Low Input indexes.
  - In column 7 under **I5\_Index\_ID**, enter the name of the SureSelect XT Low Input Dual Index P5 Indexed Adaptor assigned to the sample. In column 8 under **index2**, enter the corresponding P5 index sequence. See [Table 53](#) on page 99 for nucleotide sequences of the index segment of the SureSelect XT Low Input Dual Index P5 Indexed Adaptors.

[Header]									
IEMFileVe	5								
Experim	XT_Low_Input								
Date	#####								
Workflow	GenerateFASTQ								
Applicatio	FASTQ Only								
Instrumer	MiSeq								
Assay	TruSeq Nano DNA								
Index Ada	TruSeq DNA CD Indexes (96 Indexes)								
Description									
Chemistry	Amplicon								
[Reads]									
	100								
	100								
[Settings]									
ReverseCo	0								
[Data]									
Sample_ID	Sample_N	Sample_P	Sample_V	Index_Pla	I7_Index_ID	index	I5_Index_ID	index2	Sample
Sample_1	Sample1	Plate1	A01	A01	A01	GTCTGTCA	A01	CAACGAGC	
Sample_2	Sample2	Plate1	B01	B01	B01	TGAAGAGA	B01	GTCGACAA	
Sample_3	Sample3	Plate1	C01	C01	C01	TTCACGCA	C01	AAGAGCCT	

**Figure 15** Sample sheet for SureSelect XT Low Input dual indexed library sequencing

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

## Sequence analysis resources

Guidelines are provided below for typical NGS analysis pipeline steps appropriate for SureSelect XT Low Input dual-indexed library data analysis. Your NGS analysis pipeline may vary.

Use Illumina's bcl2fastq, BCL Convert or DRAGEN software to generate paired end reads by demultiplexing sequences based on the dual indexes. The demultiplexed FASTQ data is pre-processed to remove sequencing adaptors by each of the analysis tools described below.

### Using Agilent's Alissa Reporter software for germline DNA workflows

Alissa Reporter software provides a complete FASTQ-to-Result solution for Agilent's SureSelect assays, processing NGS data from FASTQ format to VCF format, and reporting human germline SNV, InDel and CNV calls.

Alissa Reporter is a cloud-based, multi-tenant software as a service (SaaS) product, delivering integrated pre-processing of SureSelect XT Low Input DNA library reads (adaptor trimming and deduplication) along with secondary data analysis and quality control (QC) analytics using a built-in dashboard. To obtain more information and to purchase access to the software please visit the [Alissa Reporter page at www.agilent.com](http://www.agilent.com).

Key considerations for SureSelect XT Low Input dual-indexed DNA assay steps prior to Alissa Reporter software analysis are summarized below:

- Alissa Reporter applications are available for germline analysis of human DNA libraries enriched using a pre-designed or custom SureSelect human probe (see [page 14](#)). Libraries enriched using SureSelect XT HS Human All Exon V7 or V8 probes are analyzed with the corresponding *Human All Exon V7 Germline* or *Human All Exon V8 Germline* application in Alissa Reporter. Libraries enriched using other probes, including additional pre-designed probes, are analyzed using an Alissa Reporter *Custom* application. The Alissa Reporter console provides tools for importing both pre-designed probe and custom probe designs from SureDesign and setting up a new *Custom* application for each imported design.

#### NOTE

Human All Exon V8+UTR and Human All Exon V8+NCV designs must be imported into Alissa Reporter for use as *Custom*-type applications. Use the *Catalog*-type Alissa Reporter applications, including the *Human All Exon V8 Germline* application, only for the specific probe indicated for the application without any additional design content.

## 5 Post-Capture Sample Processing for Multiplexed Sequencing

### Sequence analysis resources

- Analysis of FFPE-derived or other DNA samples for detection of somatic variants is not supported at the time of this publication. Please visit the [Alissa Reporter page at www.agilent.com](http://www.agilent.com) for information on the latest Alissa Reporter software version capabilities.
- For CNV calling a co-analysis strategy is used in which unrelated samples from the same Alissa Reporter run are used to determine the reference signal for the target sample (no specific reference sample is required). At least 3 and preferably 8 (or more) unrelated samples need to be analyzed in Alissa Reporter together to obtain a reliable reference signal for CNV calls. For CNV calling on the X and Y chromosomes, unrelated samples of the same gender are required. For best results, process the samples to be used for CNV co-analysis in the same SureSelect run and in the same sequencing run in order to minimize any processing-based variance.
- Maximum file size for uploads is 50GB/file (in total 400GB/sample). A maximum of 768 FASTQ files can be uploaded in a run.
- File sizes >150M reads are randomly subsampled to 150M reads when using the *Human All Exon V7 Germline* or *Human All Exon V8 Germline* application.
- Unmerged and merged FASTQ files are supported. Upload of BAM files or other non-FASTQ file formats is not supported.
- During Alissa Reporter run setup, in the software's **Application chemistry** menu select *XTHS (no MBC)* for analysis of the dual-indexed SureSelect XT Low Input libraries.

### Using Agilent's AGeNT software for germline or somatic DNA workflows

Agilent's AGeNT software is a Java-based toolkit used for SureSelect XT Low Input DNA library read processing steps. The AGeNT tools are designed to enable building, integrating, maintaining, and troubleshooting internal analysis pipelines for users with bioinformatics expertise. For additional information and to download this toolkit, visit the [AGeNT page at www.agilent.com](http://www.agilent.com) and review [AGeNT Best Practices](#) for processing steps suitable for SureSelect XT HS/Low Input DNA libraries, summarized below.

Prior to variant discovery, demultiplexed FASTQ data are pre-processed using the AGeNT Trimmer module to remove sequencing adaptors. The trimmed reads should be aligned using a suitable tool such as BWA-MEM. Once alignment is complete, the AGeNT CReaK (Consensus Read Kit) tool

is used to generate consensus reads and mark or remove duplicates. The resulting BAM files are ready for downstream analysis including variant discovery.

**NOTE**

CReaK is a deduplication tool introduced in AGeNT version 3.0, replacing the AGeNT LocatIt tool. Please visit the [AGeNT page at www.agilent.com](http://www.agilent.com) and review the FAQs for a detailed comparison of LocatIt and CReaK. LocatIt remains available for backward compatibility but CReaK is the recommended tool.

---

**5 Post-Capture Sample Processing for Multiplexed Sequencing**  
Sequence analysis resources



## 6 Appendix: Using FFPE-derived DNA Samples

Protocol modifications for FFPE Samples	88
Methods for FFPE Sample Qualification	88
Sequencing Output Recommendations for FFPE Samples	89

This chapter summarizes the protocol modifications to apply to FFPE samples based on the integrity of the FFPE sample DNA.



## Protocol modifications for FFPE Samples

Protocol modifications that should be applied to FFPE samples are summarized in [Table 39](#).

**Table 39** Summary of protocol modifications for FFPE samples

Workflow Step and page	Parameter	Condition for non-FFPE Samples	Condition for FFPE Samples
gDNA Sample Preparation <a href="#">page 23</a>	Qualification of DNA Integrity	Not required	Required
DNA input for Library Preparation <a href="#">page 23</a>	Input amount and means of quantification	10 ng to 200 ng, quantified by Qubit assay	Based on determined DNA integrity (see <a href="#">Table 8</a> on page 24 and <a href="#">Table 9</a> on page 25)
DNA Shearing <a href="#">page 26</a>	Mode of DNA Shearing	2 × 120 seconds	240 seconds (continuous)
Enzymatic Fragmentation of DNA <a href="#">page 29</a>	Duration of 37°C incubation	2 × 100 reads: 15 minutes 2 × 150 reads: 10 minutes	2 × 100 reads: 15 minutes 2 × 150 reads: 15 minutes
Pre-capture PCR <a href="#">page 43</a>	Cycle number	8–11	11–14
Sequencing <a href="#">page 89</a>	Output augmentation	Per project requirements	1× to 10× based on determined DNA integrity (see <a href="#">Table 40</a> and <a href="#">Table 41</a> on page 89)

## Methods for FFPE Sample Qualification

DNA integrity may be assessed using the Agilent NGS FFPE QC Kit or using the Agilent TapeStation instrument and Genomic DNA ScreenTape.

The Agilent NGS FFPE QC Kit provides a qPCR-based assay for DNA sample integrity determination. Results include the precise quantity of amplifiable DNA in the sample to allow direct normalization of input DNA amount and a  $\Delta\Delta Cq$  DNA integrity score used to design other protocol modifications.

The Agilent TapeStation instrument, combined with the Genomic DNA ScreenTape assay, provides an automated electrophoresis method for determination of a DNA Integrity Number (DIN) score used to estimate amount of input DNA required for sample normalization and to design other protocol modifications.



## Sequencing Output Recommendations for FFPE Samples

After determining the amount of sequencing output required for intact DNA samples to meet the goals of your project, use the guidelines below to determine the amount of extra sequencing output required for FFPE DNA samples.

**Samples qualified using  $\Delta\Delta Cq$ :** For samples qualified based on the  $\Delta\Delta Cq$  DNA integrity score, use the guidelines in [Table 40](#). For example, if your workflow demands 100 Mb output for intact DNA samples to achieve the required coverage, an FFPE sample with  $\Delta\Delta Cq$  score of 1 requires 200–400 Mb of sequencing output to achieve the same coverage.

**Table 40** Recommended sequencing augmentation for FFPE-derived DNA samples

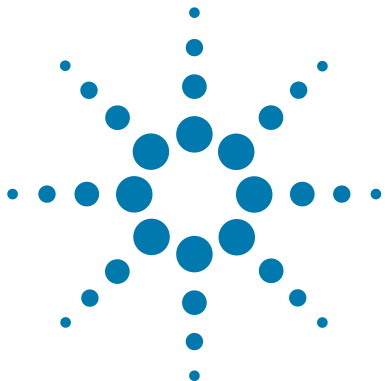
$\Delta\Delta Cq$ value	Recommended fold increase for FFPE-derived sample
<0.5	No extra sequencing output
between 0.5 and 2	Increase sequencing allocation by 2× to 4×
>2	Increase sequencing allocation by 5× to 10× or more

**Samples qualified using DIN:** For samples qualified based on the Genomic DNA ScreenTape assay DIN integrity score, use the guidelines in [Table 41](#). For example, if your workflow demands 100 Mb output for intact DNA samples to achieve the required coverage, an FFPE sample with DIN score of 4 requires approximately 200–400 Mb of sequencing output to achieve the same coverage.

**Table 41** Recommended sequencing augmentation for FFPE-derived DNA samples

DIN value	Recommended fold increase for FFPE-derived sample
$\geq 8$	No extra sequencing output
between 3 and 8	Increase sequencing allocation by 2× to 4×
<3	Increase sequencing allocation by 5× to 10× or more

**6 Appendix: Using FFPE-derived DNA Samples**  
Sequencing Output Recommendations for FFPE Samples



## 7 Reference

Kit Contents	92
Dual Indexing Primer Information	96
Troubleshooting Guide	101
Quick Reference Protocol	106

This chapter contains reference information, including component kit contents, index sequences, troubleshooting information, and a quick-reference protocol for experienced users.



## Kit Contents

Components supplied in the Agilent kits used in this protocol are detailed below.

**Table 42** Contents of SureSelect XT Low Input Dual Index P5 Indexed Adaptors 1-96 for ILM p/n 5191-4056

Kit Component	Storage Condition	Format
SureSelect XT Low Input Dual Index P5 Indexed Adaptors 1-96 for ILM	–20°C	P5 Indexed Adaptors 1 through 96 (adaptor oligos containing 8-bp P5 index sequence), provided in green plate*

\* See [Table 50](#) on page 96 for a plate map and see [Table 52](#) on page 98 or [Table 53](#) on page 99 for index sequences.

### CAUTION

The SureSelect XT Low Input Dual Index P5 Indexed Adaptors are provided in single-use aliquots. To avoid cross-contamination of libraries, use each well for a single library preparation reaction. Do not re-use any residual volume for subsequent experiments.

**Table 43** Contents of SureSelect Enzymatic Fragmentation Kit (stored at –20°C)

Kit Component	16 Reactions (p/n 5191-4079)	96 Reactions (5191-4080)
SureSelect Fragmentation Enzyme	tube with green cap	tube with green cap
5× SureSelect Fragmentation Buffer	tube with blue cap	tube with blue cap

**Table 44** Contents of SureSelect XT Low Input Reagent Kit p/n G9703A

Component Kit Name	Storage Condition	Component Kit p/n
SureSelect XT Low Input Index Primers 1–96 for ILM (Pre PCR)	–20°C	5190-6444 (see <a href="#">Table 45</a> )
SureSelect XT HS and XT Low Input Library Preparation Kit for ILM (Pre PCR)	–20°C	5500-0140 (see <a href="#">Table 46</a> )
SureSelect Target Enrichment Kit, ILM Hyb Module, Box 1 (Post PCR)	Room Temperature	5190-9687 (see <a href="#">Table 47</a> )
SureSelect XT HS and XT Low Input Target Enrichment Kit ILM Hyb Module, Box 2 (Post PCR)	–20°C	5190-9686 (see <a href="#">Table 48</a> )

The contents of each of the component kits listed in [Table 44](#) are described in the tables below.

**Table 45** SureSelect XT Low Input Index Primers for ILM Kits (Pre PCR) Content

Kit Component	Format
SureSelect XT Low Input Index Primers 1–96 for ILM	Reverse PCR primers containing 8-bp P7 index sequence 1 through 96, provided in yellow plate (Index Plate 1)*

\* See [Table 50](#) on page 96 for a plate map and see [Table 52](#) on page 98 for index sequences.

### CAUTION

The SureSelect XT Low Input Index Primers are provided in single-use aliquots. To avoid cross-contamination of libraries, use each well for a single library preparation reaction. Do not re-use any residual volume for subsequent experiments.

**Table 46** SureSelect XT HS and XT Low Input Library Preparation Kit (Pre PCR) Content

Kit Component	Format
End Repair-A Tailing Enzyme Mix	tube with orange cap
End Repair-A Tailing Buffer	bottle
T4 DNA Ligase	tube with blue cap
Ligation Buffer	bottle
Adaptor Oligo Mix*	tube with white cap
Forward Primer	tube with brown cap
100 mM dNTP Mix (25 mM each dNTP)	tube with green cap
Herculase II Fusion DNA Polymerase	tube with red cap
5× Herculase II Reaction Buffer	tube with clear cap

\* The Adaptor Oligo Mix is not used in the dual indexed library preparation protocol described in this publication.

**Table 47** SureSelect Target Enrichment Kit, ILM Hyb Module Box 1 (Post PCR) Content

Kit Component	Format
SureSelect Binding Buffer	bottle
SureSelect Wash Buffer 1	bottle
SureSelect Wash Buffer 2	bottle

**Table 48** SureSelect XT HS and XT Low Input Target Enrichment Kit, ILM Hyb Module Box 2 (Post PCR) Content

Kit Component	Format
SureSelect Fast Hybridization Buffer	bottle
SureSelect XT HS and XT Low Input Blocker Mix	tube with blue cap
SureSelect RNase Block	tube with purple cap
SureSelect Post-Capture Primer Mix	tube with clear cap
100 mM dNTP Mix (25 mM each dNTP)	tube with green cap
Herculase II Fusion DNA Polymerase	tube with red cap
5× Herculase II Reaction Buffer	tube with clear cap

Bundles of a SureSelect XT Low Input Reagent Kit with certain Target Enrichment Probes are available for purchase using the Agilent part numbers listed in Table 49. The SureSelect XT Low Input Reagent Kit included in these bundles is supplied with the same component kits listed in Table 44 on page 92.

**Table 49** Supported SureSelect XT Low Input Reagent Kit + Probe Bundles

Included SureSelect (SSel) XT Low Input Probe	Bundle part number
Custom 1–499 kb*	G9707A
Custom 0.5 –2.9 Mb*	G9707B
Custom 3–5.9 Mb*	G9707C
Custom 6–11.9 Mb*	G9707D
Custom 12–24 Mb*	G9707E
ClearSeq Comp Cancer	G9707G
Clinical Research Exome V2	G9707H
Clinical Research Exome V2 Plus	G9707J
Human All Exon V6	G9707K
Human All Exon V6 Plus	G9707L
Human All Exon V6+UTRs	G9707M
Human All Exon V7	G9707N
Human All Exon V7 Plus 1	G9707P
Human All Exon V7 Plus 2	G9707Q
Cancer All-In-One Lung	G9707R
Cancer All-In-One Solid Tumor	G9707S

\* Kits that include Custom SureSelect Cancer All-In-One panels, designed using Agilent’s SureDesign application, are ordered using these bundled custom design Agilent part numbers. Custom SureSelect Cancer All-In-One panels are designated using design IDs beginning with an ‘A’ character. (Refer to the probe vial label and the associated Certificate of Analysis to view the design ID.)

## Dual Indexing Primer Information

### Index Primer Plate Map

The plate map below shows the plate well position of each index for both P7 indexes in SureSelect XT Low Input Index Primers (yellow plate) and P5 indexes in SureSelect XT Low Input Dual Index P5 Indexed Adaptors (green plate). For each individual sample prepared using the dual indexing protocol, use the same P7 and P5 index number, originating from the same well position in the two plates. Use a different P7/P5 index pair for each sample to be multiplexed in the same sequencing reaction.

**Table 50** Plate map for P7 indexes (yellow plate) and P5 indexes (green plate)

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



## Index Nucleotide Sequences

The nucleotide sequence of the index portion of each P7 indexing primer is provided in [Table 52](#) on page 98. P7 indexes are shown in forward orientation, applicable to any of the supported Illumina platforms.

Sequences of the index portion of the P5 indexed adaptors are provided in two orientations (forward and reverse complement) in [Table 53](#) on page 99 and [Table 54](#) on page 100. The P5 index sequences are provided in both orientations to facilitate use with different Illumina sequencing platforms and sequencing run setup and management tools, as summarized in [Table 51](#). Correct representation of the P5 index orientation in sample sheets or during sequencing run setup is crucial to successful demultiplexing. Refer to Illumina support documentation and resources to determine the correct P5 index orientation for your application (combination of sequencing platform and tools, e.g., Local Run Manager and Instrument Run Setup).

**Table 51** P5 index sequencing orientation by Illumina platform

P5 Index Orientation	Platform
Forward	NovaSeq 6000 with v1.0 chemistry MiSeq HiSeq 2500
Reverse Complement*	NovaSeq 6000 with v1.5 chemistry NextSeq 500/550/1000/2000 HiSeq 3000/4000 iSeq 100 MiniSeq HiSeq X

\* Some run setup and management tools used with these platforms automatically create the reverse complement sequence for the P5 index sequence entered for the run. Be sure to consult Illumina's support documentation for the combination of platform and tools used in your pipeline to determine the correct index orientation to enter during run setup.

**7 Reference**  
Index Nucleotide Sequences

**P7 Index Sequences in Forward Orientation (use with all platforms)**

**Table 52 P7 Index 1–96 (yellow 96-well plate) Forward Sequence**

Index	Well	Sequence	Index	Well	Sequence	Index	Well	Sequence	Index	Well	Sequence
1	A01	GTCTGTCA	25	A04	CCGTGAGA	49	A07	ATGCCTAA	73	A10	ACAGCAGA
2	B01	TGAAGAGA	26	B04	GACTAGTA	50	B07	ATCATTCC	74	B10	AAGAGATC
3	C01	TTCACGCA	27	C04	GATAGACA	51	C07	AACTCACC	75	C10	CAAGACTA
4	D01	AACGTGAT	28	D04	GCTCGGTA	52	D07	AACGCTTA	76	D10	AAGACGGA
5	E01	ACCACTGT	29	E04	GGTGC GAA	53	E07	CAGCGTTA	77	E10	GCCAAGAC
6	F01	ACCTCCAA	30	F04	AACAACCA	54	F07	CTCAATGA	78	F10	CTGTAGCC
7	G01	ATTGAGGA	31	G04	CGGATTGC	55	G07	AATGTTGC	79	G10	CGCTGATC
8	H01	ACACAGAA	32	H04	AGTCACTA	56	H07	CAAGGAGC	80	H10	CAACCACA
9	A02	GCGAGTAA	33	A05	AAACATCG	57	A08	GAATCTGA	81	A11	CCTCCTGA
10	B02	GTCGTAGA	34	B05	ACGTATCA	58	B08	GAGCTGAA	82	B11	TCTTCACA
11	C02	GTGTTCTA	35	C05	CCATCCTC	59	C08	GCCACATA	83	C11	GAACAGGC
12	D02	TATCAGCA	36	D05	GGAGAACA	60	D08	GCTAACGA	84	D11	ATTGGCTC
13	E02	TGGAACAA	37	E05	CGAACTTA	61	E08	GTACGCAA	85	E11	AAGGACAC
14	F02	TGGTGGTA	38	F05	ACAAGCTA	62	F08	TCCGTCTA	86	F11	ACACGACC
15	G02	ACTATGCA	39	G05	CTGAGCCA	63	G08	CAGATCTG	87	G11	ATAGCGAC
16	H02	CCTAATCC	40	H05	ACATTGGC	64	H08	AGTACAAG	88	H11	CCGAAGTA
17	A03	AGCAGGAA	41	A06	CATACCAA	65	A09	AGGCTAAC	89	A12	CCTCTATC
18	B03	AGCCATGC	42	B06	CAATGGAA	66	B09	CGACTGGA	90	B12	AACCGAGA
19	C03	TGGCTTCA	43	C06	ACGCTCGA	67	C09	CACCTTAC	91	C12	GATGAATC
20	D03	CATCAAGT	44	D06	CCAGTTCA	68	D09	CACTTCGA	92	D12	GACAGTGC
21	E03	CTAAGGTC	45	E06	TAGGATGA	69	E09	GAGTTAGC	93	E12	CCGACAAC
22	F03	AGTGGTCA	46	F06	CGCATACA	70	F09	CTGGCATA	94	F12	AGCACCTC
23	G03	AGATCGCA	47	G06	AGAGTCAA	71	G09	AAGGTACA	95	G12	ACAGATTC
24	H03	ATCCTGTA	48	H06	AGATGTAC	72	H09	CGACACAC	96	H12	AATCCGTC

**P5 Index Sequences in Forward Orientation (see Table 51 for usage guidelines)**

**Table 53 P5 Index 1–96 (green 96-well plate) Forward Sequence**

Index	Well	Sequence	Index	Well	Sequence	Index	Well	Sequence	Index	Well	Sequence
1	A01	CAACGAGC	25	A04	CCTGCGTG	49	A07	CGATACAG	73	A10	ACTCAGGC
2	B01	GTCGACAA	26	B04	TCCAACA	50	B07	CCGTA CTC	74	B10	GACCGCAT
3	C01	AAGAGCCT	27	C04	AGGCCTAG	51	C07	GTCGGTGT	75	C10	AGAGAGAA
4	D01	ACACCTTA	28	D04	AACGTGTC	52	D07	CTGACTAA	76	D10	TCAGATTG
5	E01	TGATCGCG	29	E04	GAGCCGCT	53	E07	ACTGGATG	77	E10	AAGATAGC
6	F01	TGGCTAGA	30	F04	ACATTACG	54	F07	TTGGCATG	78	F10	CGATTGGT
7	G01	GCCTCCGA	31	G04	TGCGACAT	55	G07	GACTCGTT	79	G10	GTTCCAAG
8	H01	CAGCGTTG	32	H04	TGTCCGGC	56	H07	CTCCGAAC	80	H10	CAATCTCG
9	A02	GTGTCTCA	33	A05	CTGTTCGC	57	A08	AATGGCAT	81	A11	CAAGTCAA
10	B02	ATAACATC	34	B05	TCACGCGA	58	B08	TAGCTGTA	82	B11	CAGTCGTG
11	C02	AACTTCCT	35	C05	GTTGTTCT	59	C08	GCTCACAC	83	C11	TTACAGTG
12	D02	GCGTTGGT	36	D05	ATTAACCG	60	D08	CTAATGTT	84	D11	ACCGGCCT
13	E02	CTAGCAAC	37	E05	CGTAGTAA	61	E08	CGTTACGT	85	E11	TCATTCCA
14	F02	TCTCGATC	38	F05	CACGCTGT	62	F08	GCGCATCA	86	F11	GCTAGGAT
15	G02	GTATGCGC	39	G05	TGGAACAG	63	G08	CCATCTAA	87	G11	ATGAATTG
16	H02	AGGTCGTT	40	H05	GTGTCCGC	64	H08	CGTCTCTT	88	H11	TTAGGCTC
17	A03	GTC AATAG	41	A06	AGAGTTCG	65	A09	ACCTTGTT	89	A12	TAACACCA
18	B03	CCTGTGAC	42	B06	TAGAACGC	66	B09	TATCGACG	90	B12	ACACTCTT
19	C03	GAGGAATA	43	C06	GAGATTAT	67	C09	TTGGCGAC	91	C12	CTGTATGA
20	D03	TGCTATCT	44	D06	TAATGAGA	68	D09	CGGAAGAT	92	D12	TTGGTCAA
21	E03	GATATCAC	45	E06	CTTGCCAA	69	E09	CAAGTATT	93	E12	CGTTGGCA
22	F03	CCTGAAGA	46	F06	CGCACAGA	70	F09	TGACGACT	94	F12	TCCACTTG
23	G03	TCTCTCAA	47	G06	GCGACTGT	71	G09	CGGCCATA	95	G12	AACGGTCA
24	H03	TTCCGTCT	48	H06	AGAATAAC	72	H09	TAAGTGGT	96	H12	CTGGACCA

**7 Reference**  
Index Nucleotide Sequences

**P5 Index Sequences in Reverse Orientation (see Table 51 for usage guidelines)**

**Table 54 P5 Index 1–96 (green 96-well plate) Reverse Sequence**

Index	Well	Sequence	Index	Well	Sequence	Index	Well	Sequence	Index	Well	Sequence
1	A01	GCTCGTTG	25	A04	CACGCAGG	49	A07	CTGTATCG	73	A10	GCCTGAGT
2	B01	TTGTGAC	26	B04	TGTTGGAA	50	B07	GAGTACGG	74	B10	ATGCGGTC
3	C01	AGGCTCTT	27	C04	CTAGGCCT	51	C07	ACACCGAC	75	C10	TTCTCTCT
4	D01	TAAGGTGT	28	D04	GACACGTT	52	D07	TTAGTCAG	76	D10	CAATCTGA
5	E01	CGCGATCA	29	E04	AGCGGCTC	53	E07	CATCCAGT	77	E10	GCTATCTT
6	F01	TCTAGCCA	30	F04	CGTAATGT	54	F07	CATGCCAA	78	F10	ACCAATCG
7	G01	TCGGAGGC	31	G04	ATGTCGCA	55	G07	AACGAGTC	79	G10	CTTGGAAC
8	H01	CAACGCTG	32	H04	GCCGGACA	56	H07	GTTGCGAG	80	H10	CGAGATTG
9	A02	TGAGACAC	33	A05	GCGAACAG	57	A08	ATGCCATT	81	A11	TTGACTTG
10	B02	GATGTTAT	34	B05	TCGCGTGA	58	B08	TACAGCTA	82	B11	CACGACTG
11	C02	AGGAAGTT	35	C05	AGAACAAC	59	C08	GTGTGAGC	83	C11	CACTGTAA
12	D02	ACCAACGC	36	D05	CGGTTAAT	60	D08	AACATTAG	84	D11	AGGCCGGT
13	E02	GTTGCTAG	37	E05	TTACTACG	61	E08	ACGTAACG	85	E11	TGGAATGA
14	F02	GATCGAGA	38	F05	ACAGCGTG	62	F08	TGATGCGC	86	F11	ATCCTAGC
15	G02	GCGCATAAC	39	G05	CTGTTCCA	63	G08	TTAGATGG	87	G11	CAATTCAT
16	H02	AACGACCT	40	H05	GCCGACAC	64	H08	AAGAGACG	88	H11	GAGCCTAA
17	A03	CTATTGAC	41	A06	CGAACTCT	65	A09	AACAAGGT	89	A12	TGGTGTTA
18	B03	GTCACAGG	42	B06	GCGTTCTA	66	B09	CGTCGATA	90	B12	AAGAGTGT
19	C03	TATTCCTC	43	C06	ATAATCTC	67	C09	GTCGCCAA	91	C12	TCATACAG
20	D03	AGATAGCA	44	D06	TCTCATTA	68	D09	ATCTCCG	92	D12	TTGACCAA
21	E03	GTGATATC	45	E06	TTGGCAAG	69	E09	AATACTTG	93	E12	TGCCAACG
22	F03	TCTTCAGG	46	F06	TCTGTGCG	70	F09	AGTCGTCA	94	F12	CAAGTGGA
23	G03	TTGAGAGA	47	G06	ACAGTCGC	71	G09	TATGGCCG	95	G12	TGACCGTT
24	H03	AGACGGAA	48	H06	GTTATTCT	72	H09	ACCACTTA	96	H12	TGGTCCAG

## Troubleshooting Guide

### If recovery of gDNA from samples is low

- ✓ Using excess tissue for gDNA isolation can reduce yield. Use only the amount of each specific tissue type recommended by the gDNA isolation protocol.
- ✓ Tissue sample lysis may not have been optimal during gDNA isolation. Monitor the extent of sample lysis during the Proteinase K digestion at 56°C by gently pipetting the digestion reaction every 20–30 minutes, visually inspecting the solution for the presence of tissue clumps. If clumps are still present after the 1-hour incubation at 56°C, add another 10 µl of Proteinase K and continue incubating at 56°C, with periodic mixing and visual inspections, for up to two additional hours. When the sample no longer contains clumps of tissue, move the sample to room temperature until lysis is complete for the remaining samples. Do not over-digest. Individual samples may be kept at room temperature for up to 2 hours before resuming the protocol. Do not exceed 3 hours incubation at 56°C for any sample.

### If concentration of FFPE DNA samples is too low for enzymatic fragmentation

- ✓ The standard enzymatic fragmentation protocol requires 10–200 ng input DNA in a volume of 7 µl, and uses a final fragmentation reaction volume of 10 µl. For dilute FFPE samples, enzymatic fragmentation may be performed using the modified protocol below:
  - Bring FFPE samples containing 10–200 ng DNA to 17 µl final volume with 1X Low TE Buffer.
  - Prepare the Fragmentation master mix as directed in [Table 13](#) on page 30.
  - Add 3 µl of the master mix to each 17-µl DNA sample. Mix and spin as directed on [page 30](#).
  - Run the thermal cycling program in [Table 11](#) on page 29 using the 37°C fragmentation duration shown in the table below.

NGS read length	High-quality DNA samples	FFPE DNA samples
2 × 100 reads	25 minutes	25 minutes
2 × 150 reads	15 minutes	25 minutes

### **If yield of pre-capture libraries is low**

- ✓ The library preparation protocol includes specific thawing, temperature control, pipetting, and mixing instructions which are required for optimal performance of the highly viscous buffer and enzyme solutions used in the protocol. Be sure to adhere to all instructions when setting up the reactions.
- ✓ Ensure that the ligation master mix (see [page 35](#)) is kept at room temperature for 30–45 minutes before use.
- ✓ PCR cycle number may require optimization. Repeat library preparation for the sample, increasing the pre-capture PCR cycle number by 1 to 2 cycles. If a high molecular weight peak (>500 bp) is observed in the electropherogram for a sample with low yield, the DNA may be overamplified. Repeat library preparation for the sample, decreasing the pre-capture PCR cycle number by 1 to 3 cycles.
- ✓ DNA isolated from degraded samples, including FFPE tissue samples, may be over-fragmented or have modifications that adversely affect library preparation processes. Use the Agilent NGS FFPE QC Kit to determine the precise quantity of amplifiable DNA in the sample and allow direct normalization of input DNA amount.
- ✓ Performance of the solid-phase reversible immobilization (SPRI) purification step may be poor. Verify the expiration date for the vial of AMPure XP beads used for purification. Adhere to all bead storage and handling conditions recommended by the manufacturer. Ensure that the beads are kept at room temperature for at least 30 minutes before use. Use freshly-prepared 70% ethanol for each SPRI procedure.
- ✓ DNA elution during SPRI purification steps may be incomplete. Ensure that the AMPure XP beads are not overdried just prior to sample elution.

### **If solids observed in the End Repair-A Tailing Buffer**

- ✓ Vortex the solution at high speed until the solids are dissolved. The observation of solids when first thawed does not impact performance, but it is important to mix the buffer until all solutes are dissolved.

### **If pre-capture library fragment size is larger than expected in electropherograms**

- ✓ Shearing may not be optimal. For intact, high-quality DNA samples, ensure that shearing is completed using the two-round shearing protocol provided, including all spinning and vortexing steps.

- ✓ Any bubbles present on the microTUBE filament may disrupt complete shearing. Spin the microTUBE for 30 seconds before the first round of shearing to ensure that any bubbles are released.

### **If pre-capture library fragment size is different than expected in electropherograms**

- ✓ FFPE DNA pre-capture libraries may have a smaller fragment size distribution due to the presence of DNA fragments in the input DNA that are smaller than the target DNA shear size.
- ✓ DNA fragment size selection during SPRI purification depends upon using the correct ratio of sample to AMPure XP beads. Before removing an aliquot of beads for the purification step, mix the beads until the suspension appears homogeneous and consistent in color and verify that you are using the bead volume recommended for pre-capture purification on [page 45](#).

### **If low molecular weight adaptor-dimer peak is present in pre-capture library electropherograms**

- ✓ The presence of a low molecular weight peak, in addition to the expected peak, indicates the presence of adaptor-dimers in the library. It is acceptable to proceed to target enrichment with library samples for which adaptor-dimers are observed in the electropherogram at low abundance, similar to the samples analyzed on [page 48](#) to [page 51](#). The presence of excessive adaptor-dimers in the samples may be associated with reduced yield of pre-capture libraries. If excessive adaptor-dimers are observed, verify that the adaptor ligation protocol is being performed as directed on [page 39](#). In particular, ensure that the Ligation master mix is mixed with the sample prior to adding the SureSelect XT Low Input Dual Index P5 Indexed Adaptor to the mixture. Do not add the Ligation master mix and the P5 Indexed Adaptor to the sample in a single step.
- ✓ For whole-genome sequencing (not specifically supported by this protocol), samples with an adaptor-dimer peak must be subjected to an additional round of SPRI-purification. To complete, first dilute the sample to 50 µl with nuclease free water, then follow the SPRI purification procedure on [page 45](#).

### If yield of post-capture libraries is low

- ✓ PCR cycle number may require optimization. Repeat library preparation and target enrichment for the sample, increasing the post-capture PCR cycle number by 1 to 2 cycles.
- ✓ The RNA Probe used for hybridization may have been compromised. Verify the expiration date on the Probe vial or Certificate of Analysis. Adhere to the recommended storage and handling conditions. Ensure that the Probe Hybridization Mix is prepared immediately before use, as directed on [page 57](#), and that solutions containing the Probe are not held at room temperature for extended periods.

### If post-capture library fragment size is different than expected in electropherograms

- ✓ DNA fragment size selection during SPRI purification depends upon using the correct ratio of sample to AMPure XP beads. Before removing an aliquot of beads for the purification step, mix the beads until the suspension appears homogeneous and consistent in color and verify that you are using the bead volume recommended for post-capture purification on [page 67](#).

### If low % on-target is observed in library sequencing results

- ✓ Stringency of post-hybridization washes may have been lower than required. Complete the wash steps as directed, paying special attention to the details of SureSelect Wash Buffer 2 washes listed below:
  - SureSelect Wash Buffer 2 is pre-warmed to 70°C (see [page 60](#)). Select a thermal cycler with a block configured for efficient heating of 0.2 ml liquid volumes; ensure that the plasticware containing the wash buffer is fully seated in the thermal cycler block wells, with minimal liquid volume visible above the block during the pre-warming step.
  - Samples are maintained at 70°C during washes (see [page 61](#))
  - Bead suspensions are mixed thoroughly during washes by pipetting up and down **and** vortexing (see [page 61](#))
- ✓ Minimize the amount of time that hybridization reactions are exposed to RT conditions during hybridization setup. Locate a vortex and plate spinner or centrifuge in close proximity to thermal cycler to retain the 65°C sample temperature during mixing and transfer steps ([step 8](#) to [step 9](#) on [page 58](#)).



**If low uniformity of coverage with high AT-dropout is observed in library sequencing results**

- ✓ High AT-dropout may indicate that hybridization conditions are too stringent to obtain the desired level of coverage for AT-rich targets. Repeat target enrichment at lower stringency using a modified thermal cycler program for hybridization, reducing the hybridization temperature in segments 4 and 5 from 65°C to 62.5°C or 60°C (see [Table 26](#) on page 55).

## Quick Reference Protocol

An abbreviated summary of the protocol steps is provided below for experienced users. Use the complete protocol on [page 22](#) to [page 82](#) until you are familiar with all of the protocol details such as reagent mixing instructions and instrument settings.

Step	Summary of Conditions
<b>Library Prep</b>	
Prepare, qualify, and fragment DNA samples	Prepare 10–200 ng gDNA (in 50 µl Low TE for Covaris or in 7 µl H <sub>2</sub> O for enzymatic fragmentation) For FFPE DNA, qualify integrity and adjust input amount as directed on <a href="#">page 24</a> and <a href="#">page 25</a> Mechanically shear DNA using Covaris with shearing conditions on <a href="#">page 27</a> OR fragment DNA using SureSelect Enzymatic Fragmentation Kit with protocol on <a href="#">page 29</a> (50 µl final volume)
Prepare Ligation master mix	Per reaction: 23 µl Ligation Buffer + 2 µl T4 DNA Ligase Keep at room temperature 30–45 min before use
Prepare End-Repair/dA-Tailing master mix	Per reaction: 16 µl End Repair-A Tailing Buffer + 4 µl End Repair-A Tailing Enzyme Mix Keep on ice
End-Repair and dA-Tail the sheared DNA	50 µl fragmented DNA sample + 20 µl End Repair/dA-Tailing master mix Incubate in thermal cycler: 15 min @ 20°C, 15 min @ 72°C, Hold @ 4°C
Ligate P5 indexed adaptor	70 µl DNA sample + 25 µl Ligation master mix + 5 µl assigned P5 Indexed Adaptor (green plate) Incubate in thermal cycler: 30 min @ 20°C, Hold @ 4°C
Purify DNA	100 µl DNA sample + 80 µl AMPure XP bead suspension Elute DNA in 35 µl nuclease-free H <sub>2</sub> O
Prepare PCR master mix	Per reaction: 10 µl 5× Herculase II Reaction Buffer + 0.5 µl 100 mM dNTP Mix + 2 µl Forward Primer + 1 µl Herculase II Fusion DNA Polymerase Keep on ice
Amplify the purified DNA and add P7 index	34.5 µl purified DNA + 13.5 µl PCR master mix + 2 µl assigned SureSelect XT Low Input Index Primer (yellow plate) Amplify in thermal cycler using program on <a href="#">page 43</a>
Purify amplified DNA	50 µl amplified DNA + 50 µl AMPure XP bead suspension Elute DNA in 15 µl nuclease-free H <sub>2</sub> O
Quantify and qualify DNA	Analyze using Agilent 2100 Bioanalyzer or 4200/4150 TapeStation instrument

Step	Summary of Conditions
<b>Hybridization/Capture</b>	
Program thermal cycler	Input thermal cycler program on <a href="#">page 55</a> and pause program
Prep DNA in hyb plate	Adjust 500–1000 ng purified prepared library to 12 µl volume with nuclease-free H <sub>2</sub> O
Run pre-hybridization blocking protocol	12 µl library DNA + 5 µl SureSelect XT HS and XT Low Input Blocker Mix Run paused thermal cycler program segments 1 through 3; start new pause during segment 3 (1 min @ 65°C)
Prepare Probe Hyb Mix	Prepare 25% RNase Block dilution, then prepare appropriate mixture below: <b>Probes ≥3 Mb:</b> 2 µl 25% RNase Block + 5 µl Probe + 6 µl SureSelect Fast Hybridization Buffer <b>Probes &lt;3 Mb:</b> 2 µl 25% RNase Block + 2 µl Probe + 3 µl nuclease-free H <sub>2</sub> O + 6 µl SureSelect Fast Hybridization Buffer
Run the hybridization	With cycler paused and samples retained in cycler, add 13 µl Probe Hyb Mix to wells Resume the thermal cycler program, completing the remaining hybridization segment and 65°C or 21°C hold segment
Prepare streptavidin beads	Wash 50 µl Dynabeads MyOne Streptavidin T1 beads 3× in 200 µl SureSelect Binding Buffer
Capture hybridized libraries	Add hybridized samples (~30 µl) to washed streptavidin beads (200 µl) Incubate 30 min at RT with vigorous shaking (1400-1800 rpm) During incubation, pre-warm 6 × 200 µl aliquots per sample of SureSelect Wash Buffer 2 to 70°C
Wash captured libraries	Collect streptavidin beads with magnetic stand, discard supernatant Wash beads 1× with 200 µl SureSelect Wash Buffer 1 at RT Wash beads 6× with 200 µl pre-warmed SureSelect Wash Buffer 2 (5 minutes at 70°C per wash) Resuspend washed beads in 25 µl nuclease-free H <sub>2</sub> O
<b>Post-capture amplification</b>	
Prepare PCR master mix	Per reaction: 12.5 µl nuclease-free H <sub>2</sub> O + 10 µl 5× Herculase II Reaction Buffer + 0.5 µl 100 mM dNTP Mix + 1 µl SureSelect Post-Capture Primer Mix + 1 µl Herculase II Fusion DNA Polymerase Keep on ice
Amplify the bead-bound captured libraries	25 µl DNA bead suspension + 25 µl PCR master mix Amplify in thermal cycler using conditions on <a href="#">page 65</a>
Purify amplified DNA	Remove streptavidin beads using magnetic stand; retain supernatant 50 µl amplified DNA + 50 µl AMPure XP bead suspension Elute DNA in 25 µl nuclease-free H <sub>2</sub> O
Quantify and qualify DNA	Analyze using Agilent 2100 Bioanalyzer or 4200/4150 TapeStation instrument

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

## In This Book

This guide contains information to run the SureSelect<sup>XT</sup> Low Input target enrichment protocol with dual indexing.

© Agilent Technologies, Inc. 2018-2022

Version E0, September 2022



p/n G9703-90050



**Agilent Technologies**