



# **SureSelect<sup>QXT</sup> Whole Genome Library Prep for Illumina Multiplexed Sequencing**

**Featuring Transposase-Based  
Library Prep Technology**

## **Protocol**

Version F0, January 2021

**SureSelect platform manufactured with Agilent  
SurePrint Technology**

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



**Agilent Technologies**

## Notices

© Agilent Technologies, Inc. 2014, 2015, 2018, 2021

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

G9682-90000

### Edition

Version F0, January 2021

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. 6,627,424 and EP Pat. No.1 283 875 81, 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 describes an optimized protocol for Illumina paired-end multiplexed whole-genome library preparation using the SureSelect<sup>QXT</sup> system.

If you wish to prepare target-enriched libraries using the SureSelect<sup>QXT</sup> system, instead see publication part number G9681-90000 at [www.agilent.com](http://www.agilent.com).

### 1 Before You Begin

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

### 2 Sample Preparation

This chapter describes the steps to prepare dual-indexed gDNA sequencing libraries for the Illumina platform.

### 3 Reference

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

## What's New in Version F0

- Updates to thermal cycler and plasticware recommendations (see [Table 2](#) on page 13 and see [step 4](#) on [page 17](#))
- Updates to ordering information for AMPure XP Kits and 1X Low TE Buffer (see [Table 1](#) on page 12) and for Qubit Fluorometer (see [Table 2](#) on page 13)
- Minor updates to Agilent Bioanalyzer assay use instructions and reference document links (see [page 27](#))
- Updates to downstream sequencing support information including sequencing kit selection and seeding concentration updates (see [Table 9](#) on page 31) and support for the NovaSeq platform (see [page 31](#) through [page 36](#) and see [page 42](#))
- Updates to instructions for adaptor trimming using Agilent's AGE<sup>NT</sup> Trimmer utility and removal of SureCall adaptor trimming information (see [page 36](#) to [page 37](#))
- Removal of reference information for expired SureSelect<sup>QXT</sup> Reagent Kits p/n G9682A/G9682B, replaced by G9684A/G9684B in 2018 (see [Table 1](#) on page 12 and see [Table 21](#) through [Table 23](#) on page 40 for current Reagent Kit information)
- Updates to Technical Support contact information (see [page 2](#))
- Updates to *Notice to Purchaser* (see [page 2](#))

## What's New in Version E0

- Support for replacement of SureSelect<sup>QXT</sup> Reagent Kits p/n G9682A/G9682B with p/n G9684A/G9684B for use with Illumina's HiSeq and MiSeq platforms (see [Table 1](#) on page 12, [Table 21](#) on page 40, and [Table 19](#) on page 41 including footnotes for each table)
- Support for use of Illumina's HiSeq 3000 and HiSeq 4000 platforms for downstream sequencing steps (see [page 31](#) through [page 33](#))

- Updates to sequencing kit selection and seeding concentration guidelines (see [page 31](#))
- Updates to custom sequencing primer dilution instructions (See [page 32](#) to [page 34](#))
- Update to sequencing run setup recommendations (see “[HiSeq or NextSeq 500 platform sequencing run setup and adaptor trimming guidelines](#)” on [page 36](#))
- Updates to dual index multiplexing guidelines (see [Table 28](#) on [page 44](#))
- Updates to Agilent 2100 Bioanalyzer system ordering information (see [page 13](#))
- Updates to supplier name for materials purchased from Thermo Fisher Scientific (see [Table 1](#) on [page 12](#) and [Table 2](#) on [page 13](#))
- Addition of product guarantee and support statement (see *Note* on [page 9](#))
- Updates to Technical Support contact information (see [page 2](#))

# Content

<b>1</b>	<b>Before You Begin</b>	<b>9</b>
	Procedural Notes	10
	Safety Notes	11
	Required Reagents	12
	Required Equipment	13
<b>2</b>	<b>Sample Preparation</b>	<b>15</b>
	Step 1. Fragment and adaptor-tag the genomic DNA samples	16
	Step 2. Purify the adaptor-tagged library using AMPure XP beads	20
	Step 3. Amplify and index the adaptor-tagged DNA library	22
	Step 4. Purify the amplified library with AMPure XP beads	25
	Step 5. Assess library DNA quantity and quality using the 2100 Bioanalyzer and High Sensitivity DNA Assay	27
	Step 6. Pool samples for multiplexed sequencing	29
	Step 7. Prepare sequencing samples	31
	Step 8. Set up the sequencing run and trim adaptors from the reads	36
<b>3</b>	<b>Reference</b>	<b>39</b>
	Kit Contents	40
	Nucleotide Sequences of SureSelect <sup>OXT</sup> Dual Indexes	41
	Guidelines for Multiplexing with Dual-Indexed Samples	43

## Contents





# 1 Before You Begin

Procedural Notes	10
Safety Notes	11
Required Reagents	12
Required Equipment	13

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

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



## Procedural Notes

- The SureSelect<sup>QXT</sup> system requires high-quality DNA samples for optimal performance. Use best practices for verifying DNA sample quality before initiating the workflow. For best practice, store diluted DNA solutions at 4°C to avoid repeated freeze-thaw cycles, which may compromise DNA quality.
- Performance of the SureSelect<sup>QXT</sup> library preparation protocol is very sensitive to variations in amounts of DNA sample and other reaction components. It is important to quantify and dilute DNA samples as described on [page 17](#). Carefully measure volumes for all reaction components, and combine components as described on [page 17](#). Use best-practices for liquid handling, including regular pipette calibration, to ensure precise volume measurement.
- Use care in handling the SureSelect QXT Enzyme Mix. After removing the vial from storage at -20°C, keep on ice or in a cold block while in use. Return the vial to storage at -20°C promptly after use.
- For each protocol step that requires removal of tube cap strips, reseal the tubes with a fresh strip of domed caps. Cap deformation may result from exposure of the cap strips to the heated lid of the thermal cycler and from other procedural steps. Reuse of strip caps can cause sample loss, sample contamination, or imprecision in sample temperatures during thermal cycler incubation steps.
- 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.

- 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.
- 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.
- In general, follow Biosafety Level 1 (BSL1) safety rules.

## Safety Notes

### CAUTION

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

**1 Before You Begin**  
Required Reagents

## Required Reagents

**Table 1** Required Reagents for SureSelect<sup>OXT</sup> Library Preparation

Description	Vendor and part number
SureSelect <sup>OXT</sup> Library Prep Kit (for Illumina HiSeq, MiSeq, and NextSeq platforms)	Agilent
16 Samples	p/n G9684A
96 Samples	p/n G9684B
AMPure XP Kit	Beckman Coulter Genomics
5 mL	p/n A63880
60 mL	p/n A63881
450 mL	p/n A63882
1X Low TE Buffer (10 mM Tris-HCl, pH 7.5-8.0, 0.1 mM EDTA)	Thermo Fisher Scientific p/n 12090-015, or equivalent
100% Ethanol, molecular biology grade	Sigma-Aldrich p/n E7023
Qubit dsDNA HS Assay Kit <i>or</i>	Thermo Fisher Scientific p/n Q32851
Qubit dsDNA BR 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

## Required Equipment

**Table 2** Required Equipment for SureSelect<sup>OXT</sup> Library Preparation

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	Consult the thermal cycler manufacturer's recommendations
DNA Analysis Platform and Consumables	
Agilent 2100 Bioanalyzer Instrument	Agilent p/n G2939BA
Agilent 2100 Expert SW Laptop Bundle (optional)	Agilent p/n G2953CA
High Sensitivity DNA Kit	Agilent p/n 5067-4626
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
Centrifuge	Eppendorf Centrifuge model 5804 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
Multichannel pipette	Rainin Pipet-Lite Multi Pipette or equivalent
P10, P20, P200 and P1000 pipettes	Rainin Pipet-Lite Pipettes or equivalent
Magnetic separator*	Thermo Fisher Scientific p/n 12331D or equivalent
Vortex mixer	general laboratory supplier
Ice bucket	general laboratory supplier
Powder-free gloves	general laboratory supplier
Sterile, nuclease-free aerosol barrier pipette tips	general laboratory supplier

\* 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**  
Required Equipment



## 2 Sample Preparation

- Step 1. Fragment and adaptor-tag the genomic DNA samples 16
- Step 2. Purify the adaptor-tagged library using AMPure XP beads 20
- Step 3. Amplify and index the adaptor-tagged DNA library 22
- Step 4. Purify the amplified library with AMPure XP beads 25
- Step 5. Assess library DNA quantity and quality using the 2100 Bioanalyzer and High Sensitivity DNA Assay 27
- Step 6. Pool samples for multiplexed sequencing 29
- Step 7. Prepare sequencing samples 31
- Step 8. Set up the sequencing run and trim adaptors from the reads 36

This section contains instructions for genomic DNA sequencing library preparation for Illumina platforms.



## 2 Sample Preparation

### Step 1. Fragment and adaptor-tag the genomic DNA samples

## Step 1. Fragment and adaptor-tag the genomic DNA samples

In this step, the gDNA is enzymatically fragmented and adaptors are added to ends of the fragments in a single reaction. This step uses the SureSelect<sup>QXT</sup> Library Prep Kit components listed in [Table 3](#) in addition to some reagents obtained from other suppliers (see [Table 1](#) on page 12).

**Table 3** Reagents for DNA fragmentation and adaptor-tagging

Kit Component	Storage Location	Where Used
SureSelect QXT Stop Solution	SureSelect QXT Library Prep Kit Box 1, Room Temperature	<a href="#">page 16</a> (below)
SureSelect QXT Buffer	SureSelect QXT Library Prep Kit Box 2, –20°C	<a href="#">page 18</a>
SureSelect QXT Enzyme Mix ILM	SureSelect QXT Library Prep Kit Box 2, –20°C	<a href="#">page 18</a>

Before you begin, remove the SureSelect QXT Enzyme Mix ILM and the SureSelect QXT Buffer tubes from storage at –20°C and place on ice. Vortex each reagent vigorously to mix before use. Remove the AMPure XP beads from storage at 4°C and allow to warm up to room temperature.

#### NOTE

While obtaining components for this step, also remove the DMSO vial from the SureSelect QXT Library Prep Kit Box 2 in –20°C storage. Leave the DMSO vial at room temperature in preparation for use on [page 24](#).

For each DNA sample to be sequenced, prepare 1 library.

- 1 Verify that the SureSelect QXT Stop Solution contains 25% ethanol, by referring to the container label and the instructions below.

Before the first use of a fresh container, add 1.5 ml of ethanol to the provided bottle containing 4.5 ml of stop solution, for a final ethanol concentration of 25%. Seal the bottle then vortex well to mix. After adding the ethanol, be sure to mark the label for reference by later users.

Keep the prepared 1X SureSelect QXT Stop Solution at room temperature, tightly sealed, until it is used on [page 19](#).



## Step 1. Fragment and adaptor-tag the genomic DNA samples

- 2 Prepare reagents for the purification protocols on [page 20](#) and [page 25](#).
  - a Verify that the AMPure XP beads are being held at room temperature. The beads should be held at room temperature for at least 30 minutes before use. *Do not freeze the beads at any time.*
  - b Prepare 800  $\mu\text{l}$  of fresh 70% ethanol per sample, plus excess, for use in the purification steps. The 70% ethanol may be used for multiple steps done on the same day, when stored in a sealed container.
- 3 Quantify and dilute gDNA samples using two serial fluorometric assays:
  - a Use the Qubit dsDNA BR Assay or Qubit dsDNA HS Assay to determine the initial concentration of each gDNA sample. Follow the manufacturer's instructions for the specific assay kit and the Qubit instrument. This step is critical for successful preparation of input DNA at the required concentration to ensure optimal fragmentation.
  - b Dilute each gDNA sample with nuclease-free water to a final concentration of 50  $\text{ng}/\mu\text{l}$  in a 1.5-ml LoBind tube.
  - c Carefully measure the DNA concentration of each of the 50  $\text{ng}/\mu\text{l}$  dilutions using a second Qubit dsDNA BR or HS Assay.
  - d Adjust each gDNA sample with nuclease-free water to a final concentration of 25  $\text{ng}/\mu\text{l}$  in a 1.5-ml LoBind tube.

**CAUTION**

The duration and temperature of incubation for DNA fragmentation must be precisely controlled for optimal results. Make sure to preprogram the thermal cycler as directed in [step 4](#) before setting up the fragmentation reactions. Do not exceed 10 minutes at 45°C, as indicated in [Table 4](#).

- 4 Preprogram a thermal cycler (with the heated lid ON) with the program in [Table 4](#). Immediately pause the program, and keep paused until samples are loaded in [step 8](#).

**Table 4** Thermal cycler program for DNA fragmentation

Step	Temperature	Time
Step 1	45°C	10 minutes
Step 2	4°C	1 minute
Step 3	4°C	Hold

## 2 Sample Preparation

### Step 1. Fragment and adaptor-tag the genomic DNA samples

- 5 Before use, vortex the SureSelect QXT Buffer and SureSelect QXT Enzyme Mix ILM tubes vigorously at high speed. Note that the SSEL QXT Buffer is viscous and thorough and vigorous mixing is critical for optimal fragmentation.

These components are in liquid form when removed from  $-20^{\circ}\text{C}$  storage and should be returned to  $-20^{\circ}\text{C}$  storage promptly after use in [step 6](#).

#### CAUTION

Minor variations in volumes of the solutions combined in [step 6](#) below may result in DNA fragment size variation.

The SureSelect QXT Buffer and Enzyme Mix solutions are highly viscous. Be sure to follow the dispensing and mixing instructions in the steps below. Thorough mixing of the reagents and reactions is critical for optimal performance.

- 6 Set up the fragmentation reactions on ice using a PCR plate or strip tube. Components must be added in the order listed below. Do not pre-mix the SureSelect QXT Buffer and Enzyme Mix.
  - a To each sample well, add 17  $\mu\text{l}$  of SureSelect QXT Buffer.
  - b Add 2  $\mu\text{l}$  of each DNA sample to its assigned sample well. While dispensing the DNA, be sure to place the pipette tip at the bottom of the well.
  - c Add 1  $\mu\text{l}$  of SureSelect QXT Enzyme Mix, ILM to each sample well. While dispensing the enzyme mixture, place the pipette tip at the bottom of the well. After dispensing of the 1  $\mu\text{l}$  of enzyme mix, pipette up and down 8 to 10 times to ensure complete transfer of the viscous solution to the well.
- 7 Seal the wells, briefly spin, then mix thoroughly by vortexing the plate or strip tube at high speed for 20 seconds.
- 8 Briefly spin the samples, then immediately place the plate or strip tube in the thermal cycler and resume the thermal cycling program in [Table 4](#).
- 9 During the 10-minute incubation of samples in the thermal cycler, vigorously vortex the AMPure XP beads at high speed to ensure homogeneous distribution of beads throughout the solution so that the beads are ready for use on [page 20](#).
- 10 When the thermal cycler has completed the 1-minute incubation at  $4^{\circ}\text{C}$ , immediately place the samples on ice and proceed to [step 11](#).

## Step 1. Fragment and adaptor-tag the genomic DNA samples

**11** Add 32  $\mu$ l of 1X SureSelect QXT Stop Solution (containing 25% ethanol) to each fragmentation reaction. Seal the wells with fresh caps, then vortex at high speed for 5 seconds. Briefly spin the plate or strip tube to collect the liquid.

Incubate the samples at room temperature for 1 minute. Proceed directly to the purification protocol on [page 20](#).

## 2 Sample Preparation

### Step 2. Purify the adaptor-tagged library using AMPure XP beads

## Step 2. Purify the adaptor-tagged library using AMPure XP beads

Before you begin, verify that the AMPure XP beads have been incubated at room temperature for at least 30 minutes and that fresh 70% ethanol has been prepared for use in [step 6](#).

- 1 Verify that the AMPure XP bead suspension has been well mixed and appears homogeneous and consistent in color.
- 2 Add 52  $\mu$ l of the homogeneous bead suspension to each sample well containing the 52- $\mu$ l DNA samples. Seal the wells, then vortex for 5 seconds. Briefly spin the samples to collect the liquid, without pelleting the beads.  

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.
- 3 Incubate samples for 5 minutes at room temperature.
- 4 Put the plate or strip tube on the magnetic stand at room temperature. Wait for the solution to clear (approximately 3 to 5 minutes).
- 5 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.
- 6 Continue to keep the plate or tubes in the magnetic stand while you dispense 200  $\mu$ l of fresh 70% ethanol in each sample well.
- 7 Wait for 1 minute to allow any disturbed beads to settle, then remove the ethanol.
- 8 Repeat [step 6](#) and [step 7](#) once for a total of two washes. Make sure to remove all of the ethanol at each wash step.
- 9 Dry the samples on the thermal cycler (with lid open) at 37°C for 1 to 3 minutes. Do not overdry the samples.
- 10 Add 11  $\mu$ l of nuclease-free water to each sample well.
- 11 Seal the sample wells, then mix well on a vortex mixer and briefly spin the plate or tubes to collect the liquid.
- 12 Incubate for 2 minutes at room temperature.
- 13 Put the plate or tubes in the magnetic stand and leave for 2 minutes or until the solution in each well is clear.

## Step 2. Purify the adaptor-tagged library using AMPure XP beads

- 14 Remove each cleared supernatant (approximately 10  $\mu$ L) to wells of a fresh plate or strip tube and keep on ice. You can discard the beads at this time.

## 2 Sample Preparation

### Step 3. Amplify and index the adaptor-tagged DNA library

## Step 3. Amplify and index the adaptor-tagged DNA library

In this step, the adaptor-tagged DNA libraries are PCR amplified using an appropriate pair of dual indexing primers.

### CAUTION

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

Before you begin, thaw then vortex to mix the reagents listed in [Table 5](#). Keep all reagents except DMSO on ice.

**Table 5** Reagents for PCR amplification and indexing

Kit Component	Storage Location	Where Used
Herculase II Fusion DNA Polymerase	SureSelect QXT Library Prep Kit Box 2, -20°C	<a href="#">page 23</a>
Herculase II 5× Reaction Buffer	SureSelect QXT Library Prep Kit Box 2, -20°C	<a href="#">page 23</a>
100 mM dNTP Mix (25 mM each dNTP)	SureSelect QXT Library Prep Kit Box 2, -20°C	<a href="#">page 23</a>
SureSelect QXT P7 and P5 dual indexing primers	SureSelect QXT Library Prep Kit Box 2, -20°C	<a href="#">page 23</a>
DMSO	Transferred to Room Temperature storage on <a href="#">page 16</a>	<a href="#">page 23</a>

Prepare one indexing amplification reaction for each DNA library.

- 1 Determine the appropriate index assignments for each sample. See the [Reference](#) section for sequences of the index portion of the P7 ([page 41](#)) and P5 ([page 42](#)) indexing primers used to amplify the DNA libraries in this step.

Use a different indexing primer combination for each sample to be sequenced in the same lane.

### NOTE

For sample multiplexing, Agilent recommends maximizing index diversity on both P7 and P5 primers as required for color balance. For example, when 8-plexing, use eight different P7 index primers with two P5 index primers. See [Table 27](#) on page 43 and [Table 28](#) on page 44 for additional details.

## Step 3. Amplify and index the adaptor-tagged DNA library

- 2 Prepare the appropriate volume of PCR reaction mix, as described in [Table 6](#), on ice. Mix well on a vortex mixer and keep on ice.

**Table 6** Preparation of PCR Reaction Mix

Reagent	Volume for 1 reaction	Volume for 16 reactions (includes excess)
Nuclease-free water	22 $\mu$ L	374 $\mu$ L
Herculase II 5 $\times$ Reaction Buffer	10 $\mu$ L	170 $\mu$ L
100 mM dNTP Mix (25 mM each dNTP)	0.5 $\mu$ L	8.5 $\mu$ L
DMSO	2.5 $\mu$ L	42.5 $\mu$ L
Herculase II Fusion DNA Polymerase	1 $\mu$ L	17 $\mu$ L
<b>Total</b>	<b>36 <math>\mu</math>L</b>	<b>612 <math>\mu</math>L</b>

- 3 Add 36  $\mu$ l of the PCR reaction mixture from [step 2](#) to each 10- $\mu$ l purified library DNA sample held on ice from [step 14](#) on [page 21](#).
- 4 Add 2  $\mu$ l of the appropriate P7 dual indexing primer (P7 i1 to P7 i12) to each PCR reaction well. Add only one of the twelve possible P7 primers to each reaction well.
- 5 Add 2  $\mu$ l of the appropriate P5 dual indexing primer (P5 i13 to P5 i20) to each PCR reaction well. Add only one of the eight possible P5 primers to each reaction well.
- 6 Seal the wells and mix by vortexing gently for 5 seconds. Spin samples briefly to collect the liquid.

## 2 Sample Preparation

### Step 3. Amplify and index the adaptor-tagged DNA library

- 7 Incubate the plate or strip tube in the thermal cycler (with the heated lid ON) and run the program in [Table 7](#).

**Table 7** Thermal cycler program for PCR indexing

Segment Number	Number of Cycles	Temperature	Time
1	1	68°C	2 minutes
2	1	98°C	30 seconds
3	5	98°C	30 seconds
		56°C	30 seconds
		72°C	1 minute
4	1	4°C	Hold



## Step 4. Purify the amplified library with AMPure XP beads

Before you begin, verify that the AMPure XP beads have been kept at room temperature for at least 30 minutes and that fresh 70% ethanol has been prepared for use in [step 6](#).

- 1 Mix the AMPure XP bead suspension well so that the suspension appears homogeneous and consistent in color.
- 2 Transfer the samples to room temperature, then add 35  $\mu\text{l}$  of the homogeneous bead suspension to each sample well containing the 50- $\mu\text{l}$  amplified DNA samples. Seal the wells, then vortex for 5 seconds. Briefly spin the samples to collect the liquid.

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.

- 3 Incubate samples for 5 minutes at room temperature.
- 4 Put the plate or strip tube on the magnetic stand at room temperature. Wait for the solution to clear (approximately 3 to 5 minutes).
- 5 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.
- 6 Continue to keep the plate or tubes in the magnetic stand while you dispense 200  $\mu\text{l}$  of fresh 70% ethanol in each sample well.
- 7 Wait for 1 minute to allow any disturbed beads to settle, then remove the ethanol.
- 8 Repeat [step 6](#) and [step 7](#) once for a total of two washes. Make sure to remove all of the ethanol at each wash step.
- 9 Dry the samples on the thermal cycler (with lid open) at 37°C for 1 to 3 minutes. Do not overdry the samples.
- 10 Add 30  $\mu\text{l}$  of nuclease-free water to each sample well.
- 11 Seal the sample wells, then mix well on a vortex mixer and briefly spin the plate or tubes to collect the liquid.
- 12 Incubate for 2 minutes at room temperature.
- 13 Put the plate or tubes in the magnetic stand and leave for 2 minutes or until the solution in each well is clear.
- 14 Remove each cleared supernatant (approximately 29  $\mu\text{L}$ ) to a fresh LoBind tube. You can discard the beads at this time.

## 2 Sample Preparation

### Step 4. Purify the amplified library with AMPure XP beads

**15** Remove a 1- $\mu$ l sample of each amplified library for analysis in “Step 5. Assess library DNA quantity and quality using the 2100 Bioanalyzer and High Sensitivity DNA Assay” on page 27. Dilute each of the 1- $\mu$ l samples with 9  $\mu$ l of nuclease-free water prior to analysis.

**Stopping Point** If you do not continue to the next step, store the libraries at 4°C short term or at -20°C for long term storage.

## Step 5. Assess library DNA quantity and quality using the 2100 Bioanalyzer and High Sensitivity DNA Assay

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

### NOTE

Do not use Agilent's DNA 1000 Assay to analyze the whole genome samples. The expected distribution of whole genome library fragment sizes is not compatible with the DNA 1000 Assay. See [Figure 1](#), below, for a sample electropherogram.

The presence of magnetic beads in the samples may adversely impact the Bioanalyzer results. If you suspect bead contamination in the samples, place the plate or strip tube on the magnetic rack before withdrawing samples for analysis.

- 1 Set up the 2100 Bioanalyzer 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 a 1:10 dilution of each sample for the analysis.
- 3 Load the prepared chip into the instrument and start the run within five minutes after preparation.
- 4 Analyze the results, using the guidelines below:
  - Typical whole genome library electropherograms show a broad distribution of DNA fragments. A sample electropherogram is shown in [Figure 1](#).
  - Check the Average Size [bp] of DNA fragments reported in the Bioanalyzer results. Sequencing data may be acquired from libraries with a broad range of average fragment sizes. The protocol has been optimized, however, to produce whole genome libraries with average DNA fragment sizes between approximately 600 and 1000 bp.

### NOTE

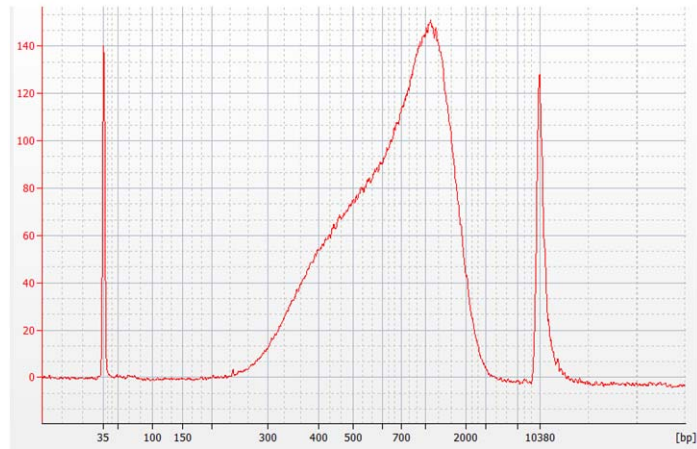
An average fragment size significantly less than 600 bp may indicate too little gDNA in the fragmentation reaction and may be associated with increased duplicates in the sequencing data. In contrast, libraries with an unusually large average fragment size may indicate too much gDNA in the fragmentation reaction and may require higher DNA concentrations for optimal cluster density in the sequencing reaction.

- 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.

## 2 Sample Preparation

### Step 5. Assess library DNA quantity and quality using the 2100 Bioanalyzer and 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.



**Figure 1** Analysis of amplified library DNA using a High-Sensitivity DNA Assay.

## Step 6. Pool samples for multiplexed sequencing

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

Guidelines for optimal low-level multiplexing of samples indexed using the SureSelect<sup>QXT</sup> dual indexes are provided on [page 43](#).

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

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

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

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

$\#$  is the number of indexes, and

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

[Table 8](#) shows an example of the amount of 4 index-tagged samples (of different concentrations) and Low TE needed for a final volume of 20  $\mu\text{L}$  at 10 nM.

**Table 8** Example of indexed sample volume calculation for total volume of 20  $\mu\text{L}$

Component	V(f)	C(i)	C(f)	#	Volume to use ( $\mu\text{L}$ )
Sample 1	20 $\mu\text{L}$	20 nM	10 nM	4	2.5
Sample 2	20 $\mu\text{L}$	10 nM	10 nM	4	5
Sample 3	20 $\mu\text{L}$	17 nM	10 nM	4	2.9
Sample 4	20 $\mu\text{L}$	25 nM	10 nM	4	2
Low TE					7.6

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

## 2 Sample Preparation

### Step 6. Pool samples for multiplexed sequencing

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

## Step 7. Prepare sequencing samples

Proceed to cluster amplification using the appropriate Illumina Paired-End Cluster Generation Kit. See [Table 9](#) for kit configurations compatible with the recommended read length plus reads for the SureSelect<sup>QXT</sup> 8-bp dual indexes.

The optimal seeding concentration for SureSelect<sup>QXT</sup> whole genome libraries varies according to sequencing platform, run type and Illumina kit version. See [Table 9](#) 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.

To do this step, refer to the manufacturer’s instructions, using the modifications described on [page 32](#) for use of the SureSelect<sup>QXT</sup> Read Primers with the Illumina Paired-End Cluster Generation Kits. Follow Illumina’s recommendation for a PhiX control in a low-concentration spike-in for improved sequencing quality control.

**Table 9** 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	14–20 pM
HiSeq 2500	High Output	2 × 100 bp	250 Cycle Kit	v4	14–20 pM
MiSeq	All Runs	2 × 100 bp or 2 × 150 bp	300 Cycle Kit	v2	14–20 pM
MiSeq	All Runs	2 × 76 bp	150 Cycle Kit	v3	14–20 pM
NextSeq 500/550	All Runs	2 × 100 bp or 2 × 150 bp	300 Cycle Kit	v2.5	2 pM
HiSeq 3000/4000	All Runs	2 × 100 bp or 2 × 150 bp	300 Cycle Kit	v1	200–300 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

\* If your application requires a different read length, verify that you have sufficient sequencing reagents to complete Reads 1 and 2 in addition to the dual 8-bp index reads.

## 2 Sample Preparation

### Step 7. Prepare sequencing samples

#### Using the SureSelect<sup>QXT</sup> Read Primers with Illumina's Paired-End Cluster Generation Kits

To sequence the SureSelect<sup>QXT</sup> libraries on Illumina's sequencing platforms, you need to use the following custom sequencing primers, provided in SureSelect QXT Library Prep Kit Box 2:

- **SureSelect QXT Read Primer 1**
- **SureSelect QXT Read Primer 2**
- **SureSelect QXT Index 1 Read Primer**
- **SureSelect QXT Index 2 Read Primer** (this primer is used only for HiSeq 3000, HiSeq 4000, and NextSeq platforms and for NovaSeq platform runs using v1.5 chemistry)

These SureSelect<sup>QXT</sup> custom sequencing primers are provided at 100  $\mu$ M and must be diluted in the corresponding Illumina primer solution, using the platform-specific instructions below:

**For the HiSeq 2500 platform**, combine the primers as shown in [Table 10](#) or [Table 11](#) on [page 33](#).

**For the HiSeq 3000 or HiSeq 4000 platform**, combine the primers as shown in [Table 12](#) on [page 33](#).

**For the MiSeq platform**, combine the primers as shown in [Table 13](#) on [page 34](#).

**For the NextSeq platform**, combine the primers as shown in [Table 14](#) or [Table 15](#) on [page 34](#).

**For the NovaSeq platform**, combine the primers as shown in [Table 16](#) through [Table 19](#) on [page 35](#).

#### NOTE

It is important to combine the primers precisely in the indicated ratios. Carefully follow the instructions indicated in [Table 10](#) to [Table 19](#). Where specified, add the custom primer volume directly to the solution already in cBot reagent plate wells. Otherwise, combine measured volumes of each solution; do not rely on volumes reported on vial labels or in Illumina literature. Vortex each mixture vigorously to ensure homogeneity for proper detection of the indexes using the custom read primers.



**Table 10 HiSeq 2500 High Output** custom sequencing primer preparation

Sequencing Read	Volume of SureSelect <sup>QXT</sup> Primer	Volume of Illumina TruSeq Primer	Total Volume
Read 1	6 µl SureSelect QXT Read Primer 1 (brown cap)	1194 µl HP10	1.2 ml*
Index	15 µl SureSelect QXT Index 1 Read Primer (clear cap)	2985 µl HP12	3 ml
Read 2	15 µl SureSelect QXT Read Primer 2 (black cap)	2985 µl HP11	3 ml

\* Aliquot the mixture as directed for HP6 or HP10 in Illumina's cluster generation protocol.

**Table 11 HiSeq 2500 Rapid Mode** custom sequencing primer preparation

Sequencing Read	Volume of SureSelect <sup>QXT</sup> Primer	Volume of Illumina TruSeq Primer	Total Volume
Read 1	8.8 µl SureSelect QXT Read Primer 1 (brown cap)	1741.2 µl HP10	1.75 ml
Index	8.8 µl SureSelect QXT Index 1 Read Primer (clear cap)	1741.2 µl HP12	1.75 ml
Read 2	8.8 µl SureSelect QXT Read Primer 2 (black cap)	1741.2 µl HP11	1.75 ml

**Table 12 HiSeq 3000 and HiSeq 4000** custom sequencing primer preparation

Sequencing Read	Volume of SureSelect <sup>QXT</sup> Primer	Volume of Illumina TruSeq Primer	Total Volume	Reagent Rack Position
Read 1	1.5 µl SureSelect QXT Read Primer 1 (brown cap)	298.5 µl HP10*	0.3 ml per well	cBot Column 11
Read 2	15 µl SureSelect QXT Read Primer 2 (black cap)	2985 µl HP11	3 ml	16
Index 1+ Index 2	22.5 µl SureSelect QXT Index 1 Read Primer (clear cap) + 22.5 µl SureSelect QXT Index 2 Read Primer (purple cap)	4455 µl HP14	4.5 ml	17

\* Use cBot recipe *HiSeq\_3000\_4000\_HD\_Exclusion\_Amp\_v1.0*. Add 1.5 µl SureSelect QXT Read Primer 1 to the 298.5 µl of HP10 in each well of column 11 in the cBot reagent plate.

## 2 Sample Preparation

### Step 7. Prepare sequencing samples

**Table 13 MiSeq platform** custom sequencing primer preparation

Sequencing Read	Volume of SureSelect <sup>QXT</sup> Primer	Volume of Illumina TruSeq Primer	Total Volume	Final Cartridge Position
Read 1	3 µl SureSelect QXT Read Primer 1 (brown cap)	597 µl HP10 (well 12)	0.6 ml	well 18
Index	3 µl SureSelect QXT Index 1 Read Primer (clear cap)	597 µl HP12 (well 13)	0.6 ml	well 19
Read 2	3 µl SureSelect QXT Read Primer 2 (black cap)	597 µl HP11 (well 14)	0.6 ml	well 20

**Table 14 NextSeq 500/550 High-Output v2 Kit** custom sequencing primer preparation

Sequencing Read	Volume of SureSelect <sup>QXT</sup> Primer	Volume of Illumina Primer	Total Volume	Final Cartridge Position
Read 1	3.9 µl SureSelect QXT Read Primer 1 (brown cap)	1296.1 µl BP10 (from well 20)	1.3 ml	well 7
Read 2	4.2 µl SureSelect QXT Read Primer 2 (black cap)	1395.8 µl BP11 (from well 21)	1.4 ml	well 8
Index 1+ Index 2	6 µl SureSelect QXT Index 1 Read Primer (clear cap) + 6 µl SureSelect QXT Index 2 Read Primer (purple cap)	1988 µl BP14 (from well 22)	2 ml	well 9

**Table 15 NextSeq 500/550 Mid-Output v2 Kit** custom sequencing primer preparation

Sequencing Read	Volume of SureSelect <sup>QXT</sup> Primer	Volume of Illumina Primer	Total Volume	Final Cartridge Position
Read 1	2.7 µl SureSelect QXT Read Primer 1 (brown cap)	897.3 µl BP10 (from well 20)	0.9 ml	well 7
Read 2	3.3 µl SureSelect QXT Read Primer 2 (black cap)	1096.7 µl BP11 (from well 21)	1.1 ml	well 8
Index 1+ Index 2	4.8 µl SureSelect QXT Index 1 Read Primer (clear cap) + 4.8 µl SureSelect QXT Index 2 Read Primer (purple cap)	1590.4 µl BP14 (from well 22)	1.6 ml	well 9

**Table 16 NovaSeq 6000 using SP/S1/S2 flowcell with v1.0 chemistry custom sequencing primer preparation**

Sequencing Read	Volume of SureSelect <sup>QXT</sup> Primer	Volume of Illumina Primer	Total Volume	Final Cartridge Position
Read 1	12 µl SureSelect QXT Read Primer 1 (brown cap)	3988 µl BP10 (well 24)	4 ml	5
Index	15 µl SureSelect QXT Index 1 Read Primer (clear cap)	4985 µl BP14 (well 23)	5 ml	7
Read 2	6 µl SureSelect QXT Read Primer 2 (black cap)	1994 µl BP11 (well 13)	2 ml	6

**Table 17 NovaSeq 6000 using S4 flowcell with v1.0 chemistry custom sequencing primer preparation**

Sequencing Read	Volume of SureSelect <sup>QXT</sup> Primer	Volume of Illumina Primer	Total Volume	Final Cartridge Position
Read 1	21.9 µl SureSelect QXT Read Primer 1 (brown cap)	7278.1 µl BP10 (well 24)	7.3 ml	5
Index	15 µl SureSelect QXT Index 1 Read Primer (clear cap)	4985 µl BP14 (well 23)	5 ml	7
Read 2	10.5 µl SureSelect QXT Read Primer 2 (black cap)	3489.5 µl BP11 (well 13)	3.5 ml	6

**Table 18 NovaSeq 6000 using SP/S1/S2 flowcell with v1.5 chemistry custom sequencing primer preparation**

Sequencing Read	Volume of SureSelect <sup>QXT</sup> Primer	Volume of Illumina Primer	Total Volume	Final Cartridge Position
Read 1	12 µl SureSelect QXT Read Primer 1 (brown cap)	3988 µl BP10 (well 24)	4 ml	5
Index 1+ Index 2	15 µl SureSelect QXT Index 1 Read Primer (clear cap) + 15 µl SureSelect QXT Index 2 Read Primer (purple cap)	4970 µl VP14 (well 23)	5 ml	7
Read 2	6 µl SureSelect QXT Read Primer 2 (black cap)	1994 µl BP11 (well 13)	2 ml	6

**Table 19 NovaSeq 6000 using S4 flowcell with v1.5 chemistry custom sequencing primer preparation**

Sequencing Read	Volume of SureSelect <sup>QXT</sup> Primer	Volume of Illumina Primer	Total Volume	Final Cartridge Position
Read 1	21.9 µl SureSelect QXT Read Primer 1 (brown cap)	7278.1 µl BP10 (well 24)	7.3 ml	5
Index 1+ Index 2	15 µl SureSelect QXT Index 1 Read Primer (clear cap) + 15 µl SureSelect QXT Index 2 Read Primer (purple cap)	4970 µl VP14 (well 23)	5 ml	7
Read 2	10.5 µl SureSelect QXT Read Primer 2 (black cap)	3489.5 µl BP11 (well 13)	3.5 ml	6

## 2 Sample Preparation

### Step 8. Set up the sequencing run and trim adaptors from the reads

## Step 8. Set up the sequencing run and trim adaptors from the reads

Refer to Illumina protocols to set up custom sequencing primer runs. Before aligning reads to the reference genome, SureSelect<sup>QXT</sup> adaptor sequences must be trimmed from the reads using the additional platform-specific guidelines below.

For SureSelect<sup>QXT</sup> dual index sequence information, see [page 41](#).

### MiSeq platform sequencing run setup and adaptor trimming guidelines

Use the Illumina Experiment Manager (IEM) software to generate a custom primer Sample Sheet.

Set up the run to include adapter trimming using the IEM Sample Sheet Wizard. When prompted by the wizard, select the *Use Adapter Trimming* option, and specify **CTGTCTCTTGATCACA** as the adaptor sequence. This enables the MiSeq Reporter software to identify the adaptor sequence and trim the adaptor from reads.

### HiSeq/NextSeq/NovaSeq platform sequencing run setup and adaptor trimming guidelines

Set up sequencing runs using the settings shown in [Table 20](#). For HiSeq runs, select *Dual Index* on the *Run Configuration* screen of the instrument control software interface. Since custom primers are spiked into the standard sequencing primer tubes, no additional specialized settings are required to accommodate the use of custom primers in the run.

For the NextSeq or NovaSeq platform, Cycle Number and custom sequencing primer settings can be specified on the *Run Configuration* screen of the instrument control software interface.

**Table 20** Run Configuration screen Cycle Number settings

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

**Step 8. Set up the sequencing run and trim adaptors from the reads**

After the sequencing run is complete, generate demultiplexed FASTQ data following Illumina's instructions and then trim adaptor sequences from the reads using the Trimmer utility of the Agilent Genomics NextGen Toolkit (AGeNT). For additional information and to download this toolkit free-of-charge, visit the [AGeNT page at www.agilent.com](http://www.agilent.com).

## 2 **Sample Preparation**

Step 8. Set up the sequencing run and trim adaptors from the reads



### 3 Reference

Kit Contents	40
Nucleotide Sequences of SureSelect <sup>OXT</sup> Dual Indexes	41
Guidelines for Multiplexing with Dual-Indexed Samples	43

This chapter contains reference information, including component kit contents and reference information for use during the downstream sample sequencing steps.



## Kit Contents

SureSelect<sup>QXT</sup> Library Prep Kits contain the following component kits:

**Table 21** SureSelect<sup>QXT</sup> Library Prep Kit Contents

Component Kits	Storage Condition	G9684A (16 Samples)	G9684B (96 Samples)
SureSelect QXT Library Prep Kit, Box 1	Room Temperature	5500-0119	5500-0119
SureSelect QXT Library Prep Kit, Box 2	-20°C	5500-0126	5500-0127

The contents of each of the component kits listed in [Table 21](#) are described in [Table 22](#) and [Table 23](#).

**Table 22** SureSelect QXT Library Prep Kit, Box 1 Content

Kit Component	16 Reactions	96 Reactions
SureSelect QXT Stop Solution	bottle	bottle

**Table 23** SureSelect QXT Library Prep Kit Box 2 Content

Kit Component	16 Reactions	96 Reactions
SureSelect QXT Buffer	tube with white cap	bottle
SureSelect QXT Enzyme Mix ILM	tube with orange cap	tube with orange cap
Herculase II Fusion DNA Polymerase	tube with red cap	tube with red cap
Herculase II 5× Reaction Buffer	tube with clear cap	tube with clear cap
100 mM dNTP Mix (25 mM each dNTP)	tube with green cap	tube with green cap
DMSO	tube with green cap	tube with green cap
SureSelect QXT Read Primer 1	tube with amber cap	tube with amber cap
SureSelect QXT Read Primer 2	tube with black cap	tube with black cap
SureSelect QXT Index 1 Read Primer	tube with clear cap	tube with clear cap
SureSelect QXT Index 2 Read Primer	tube with purple cap	tube with purple cap
SureSelect QXT P7 dual indexing primers	P7 i1 through P7 i8 provided in 8 tubes with yellow caps (one tube per primer)	P7 i1 through P7 i12 provided in 12 tubes with yellow caps (one tube per primer)
SureSelect QXT P5 dual indexing primers	P5 i13 and P5 i14 provided in 2 tubes with blue caps (one tube per primer)	P5 i13 through P5 i20 provided in 8 tubes with blue caps (one tube per primer)



## Nucleotide Sequences of SureSelect<sup>QXT</sup> Dual Indexes

The nucleotide sequence of each SureSelect<sup>QXT</sup> index is provided in the tables below.

Note that some index number assignments of the SureSelect<sup>QXT</sup> P5 and P7 indexes differ from the index numbers assignments used by Illumina for indexes of similar or identical sequence.

Each index is 8 bases in length. Refer to Illumina’s sequencing run setup instructions for sequencing libraries using 8-base indexes.

**Table 24** SureSelect<sup>QXT</sup> P7 Indexes 1 to 12

Index Name with Number	Sequence
P7 Index 1 (P7 i1)	TAAGGCCG
P7 Index 2 (P7 i2)	CGTACTAG
P7 Index 3 (P7 i3)	AGGCAGAA
P7 Index 4 (P7 i4)	TCCTGAGC
P7 Index 5 (P7 i5)	GTAGAGGA
P7 Index 6 (P7 i6)	TAGGCATG
P7 Index 7 (P7 i7)	CTCTCTAC
P7 Index 8 (P7 i8)	CAGAGAGG
P7 Index 9 (P7 i9)	GCTACGCT
P7 Index 10 (P7 i10)	CGAGGCTG
P7 Index 11 (P7 i11)	AAGAGGCA
P7 Index 12 (P7 i12)	GGACTCCT

### 3 Reference

#### Nucleotide Sequences of SureSelect<sup>OXT</sup> Dual Indexes

**Table 25** SureSelect<sup>OXT</sup> P5 Indexes 13 to 20 for HiSeq 2500, MiSeq, or NovaSeq (v1.0 chemistry) platform

Index Number	Sequence
P5 Index 13 (P5 i13)	TAGATCGC
P5 Index 14 (P5 i14)	CTCTCTAT
P5 Index 15 (P5 i15)	TATCCTCT
P5 Index 16 (P5 i16)	AGAGTAGA
P5 Index 17 (P5 i17)	GTAAGGAG
P5 Index 18 (P5 i18)	ACTGCATA
P5 Index 19 (P5 i19)	AAGGAGTA
P5 Index 20 (P5 i20)	CTAAGCCT

**Table 26** SureSelect<sup>OXT</sup> P5 Indexes 13 to 20 for HiSeq 3000/4000, NextSeq, or NovaSeq (v1.5 chemistry) platform \*

Index Number	Sequence
P5 Index 13 (P5 i13)	GCGATCTA
P5 Index 14 (P5 i14)	ATAGAGAG
P5 Index 15 (P5 i15)	AGAGGATA
P5 Index 16 (P5 i16)	TCTACTCT
P5 Index 17 (P5 i17)	CTCCTTAC
P5 Index 18 (P5 i18)	TATGCAGT
P5 Index 19 (P5 i19)	TACTCCTT
P5 Index 20 (P5 i20)	AGGCTTAG

\* When doing runs on these platforms through BaseSpace, use the reverse complement sequences provided in [Table 25](#).

## Guidelines for Multiplexing with Dual-Indexed Samples

Agilent recommends following the dual index sample pooling guidelines shown in [Table 27](#) for 16 reaction kits and shown in [Table 28](#) for 96 reaction kits. These are designed to maintain color balance at each cycle of the index reads on both ends. They also provide flexibility of demultiplexing as single or dual indexed samples in low-plexity experiments. One-base mismatches should be allowed during demultiplexing.

**Table 27** Dual index sample pooling guidelines for 16 Reaction Kits

Plexity of Sample Pool	Recommended SureSelect <sup>OXT</sup> P7 Indexes	Recommended SureSelect <sup>OXT</sup> P5 Indexes
1-plex	Any P7 index (i1 to i8)	Either P5 index (i13 or i14)
2-plex	P7 i1 and P7 i2 OR P7 i2 and P7 i4	P5 i13 and P5 i14
3-plex	P7 i1, P7 i2 and P7 i4 OR P7 i3, P7 i4 and P7 i6 OR P7 i5, P7 i7 and P7 i8	P5 i13 and P5 i14 (as needed)
4- or 5-plex	P7 i1, P7 i2, P7 i4 and any additional P7 index(es) OR P7 i3, P7 i4, P7 i6 and any additional P7 index(es) OR P7 i5, P7 i7, P7 i8 and any additional P7 index(es)	P5 i13 and P5 i14 (as needed)
6- to 8-plex	Any combination of 6, 7, or 8 different P7 indexes	P5 i13 and P5 i14 (as needed)
9-to 16-plex	All eight P7 indexes (i1 to i8)	P5 i13 and P5 i14 (as needed)

### 3 Reference

#### Guidelines for Multiplexing with Dual-Indexed Samples

**Table 28** Dual index sample pooling guidelines for 96 Reaction Kits

Plexity of Sample Pool	Recommended SureSelect <sup>QXT</sup> P7 Indexes	Recommended SureSelect <sup>QXT</sup> P5 Indexes
1-plex	Any P7 index i1 to i12	Any P5 index (i13 to i20)
2-plex	P7 i1 and P7 i2 OR P7 i2 and P7 i4	P5 i13 and P5 i14 OR P5 i15 and P5 i16 OR P5 i17 and P5 i18
3-plex	P7 i1, P7 i2 and P7 i4 OR P7 i3, P7 i4 and P7 i6 OR P7 i5, P7 i7 and P7 i8	P5 i13 and P5 i14 OR P5 i15 and P5 i16 OR P5 i17 and P5 i18 (as needed)
4-plex	P7 i1, P7 i2, P7 i3* and P7 i4 OR P7 i3, P7 i4, P7 i5* and P7 i6 OR P7 i5, P7 i6*, P7 i7 and P7 i8	P5 i13 and P5 i14 OR P5 i15 and P5 i16 OR P5 i17 and P5 i18 (as needed)
5-plex	P7 i1, P7 i2, P7 i3*, P7 i4 and P7 i5* OR P7 i3, P7 i4, P7 i5*, P7 i6 and p7 i7* OR P7 i5, P7 i6*, P7 i7, P7 i8 and p7 i9*	P5 i13 and P5 i14 OR P5 i15 and P5 i16 OR P5 i17 and P5 i18 (as needed)
6- to 12-plex	Any combination of P7 indexes i1 to i12 using each index only once	P5 i13 and P5 i14 OR P5 i15 and P5 i16 OR P5 i17 and P5 i18 (as needed)
13-to 96-plex	All twelve P7 indexes (i1 to i12)	P5 i13 and P5 i14 and any other P5 index OR P5 i15 and P5 i16 and any other P5 index OR P5 i17 and P5 i18 and any other P5 index (as needed)

\* The indicated indexes may be substituted with another index, as long as the substitute index differs from all others used in the sample pool.

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

## In This Book

This guide contains information to run the SureSelect<sup>QXT</sup> Whole Genome Library Prep protocol for Illumina paired-end multiplexed sequencing.

© Agilent Technologies, Inc. 2014, 2015,  
2018, 2021

Version F0, January 2021



p/n G9682-90000



**Agilent Technologies**