



# SureSelect Max DNA Library Preparation with Enzymatic Fragmentation

For Illumina Platform NGS

## Protocol

**Version B0 January 2026**

SureSelect platform manufactured with Agilent SurePrint technology.

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

# Notices

© Agilent Technologies, Inc. 2024, 2026

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

G9660-90000

## Edition

Version B0 January 2026

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.

## 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 Illumina paired-end multiplexed DNA sequencing libraries using the SureSelect Max Enzymatic Fragmentation Library Preparation Module. The SureSelect Max workflow segment supported by this guide includes enzymatic fragmentation of gDNA samples through library preparation using adaptors with optional duplex molecular barcodes or MBCs. Prepared libraries are ready for target enrichment as described in separate guides for later workflow segments.

### 1 Before You Begin

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

### 2 DNA Library Preparation with Enzymatic Fragmentation Protocol for Target Enrichment

This section describes the steps to prepare dual-indexed gDNA sequencing libraries using enzymatic fragmentation. Libraries can be prepared with either MBC-tagged or MBC-free adaptors. Libraries prepared using this protocol are ready for use in the SureSelect Max Target Enrichment protocols.

### 3 Appendix: Whole Genome Library Preparation Protocol

This section describes the protocol modifications required to prepare dual-indexed gDNA sequencing libraries using enzymatic fragmentation for whole genome sequencing applications. Guidelines for downstream NGS are included in this section.

### 4 Reference

This section contains reference information, including component kit contents and troubleshooting information.

## What's New in Version B0

- Added details to footnotes for [Table 1](#) on page 9 and [Table 10](#) on page 18 regarding the kit size and run configuration descriptions.
- Minor protocol updates including:
  - Revised preparative steps for the supplied reagents and user-prepared mixtures ([Table 5](#) on page 14 and instructions on [page 18](#), [page 19](#), and [page 22](#))
  - Revised *Note* on [page 19](#) regarding addition of Adaptor Oligo Mix and Ligation Master Mix
  - Clarification to FFPE DNA input amount determinants in footnotes to [Table 14](#) on page 21 and [Table 22](#) on page 29
  - Updates to guidance on solvents for sample preparation and dilution (1X Low TE Buffer or water) on [page 15](#)
- Update to the example TapeStation electropherogram in [Figure 2](#) on page 25.
- Minor updates to instructions for obtaining NGS support for SureSelect Max for ILM library sequencing guidance ([page 36](#)).
- Fully updated "[Analysis pipeline guidelines](#)" on page 36.
- New *Note* below [Table 1](#) on page 9 explaining use of AMPure XP Beads as a replacement for SureSelect Max Purification Beads.

# Content

|          |  |           |
|----------|--|-----------|
| <b>1</b> | <b>Before You Begin</b>  | <b>7</b>  |
|          | Overview of the Workflow   | <b>8</b>  |
|          | SureSelect Max Modules Used in the Workflow  | <b>9</b>  |
|          | Additional Materials Used in the Workflow  | <b>10</b> |
|          | Procedural and Safety Notes  | <b>12</b> |
| <b>2</b> | <b>DNA Library Preparation with Enzymatic Fragmentation Protocol for Target Enrichment</b> | <b>13</b> |
|          | Step 1. Prepare and qualify genomic DNA samples  | <b>15</b> |
|          | Preparation of high-quality gDNA from fresh biological samples                             | 15        |
|          | Preparation and qualification of gDNA from FFPE samples                                    | 15        |
|          | Step 2. Fragment, end-repair, and 3'-dA-tail the DNA (Frag/A-Tail)                         | <b>17</b> |
|          | Step 3. Ligate the adaptor   | <b>19</b> |
|          | Step 4. Purify libraries using magnetic purification beads                                 | <b>19</b> |
|          | Step 5. Amplify and index the libraries  | <b>21</b> |
|          | Step 6. Purify amplified libraries using magnetic purification beads                       | <b>22</b> |
|          | Step 7. QC and quantify the libraries (optional)   | <b>24</b> |
| <b>3</b> | <b>Appendix: Whole Genome Library Preparation Protocol</b>                                 | <b>27</b> |
|          | Protocol Overview  | <b>28</b> |
|          | Step 1. Prepare and qualify genomic DNA samples  | 28        |
|          | Step 2. Fragment, end-repair, and 3'-dA-tail the DNA (Frag/A-Tail)                         | 28        |
|          | Step 3. Ligate the adaptor   | 29        |
|          | Step 4. Purify libraries using magnetic purification beads                                 | 29        |
|          | Step 5. Index and amplify the libraries  | 29        |
|          | Step 6. Purify amplified libraries using magnetic purification beads                       | 30        |
|          | Step 7. QC and quantify the libraries  | 32        |
|          | NGS Guidelines   | <b>34</b> |
|          | Library pooling guidelines for multiplex NGS   | 34        |
|          | Sequencing setup and run guidelines  | 35        |
|          | Analysis pipeline guidelines   | 36        |
| <b>4</b> | <b>Reference</b>   | <b>38</b> |
|          | Reagent Kit Contents   | <b>39</b> |
|          | Component Kit Details  | 39        |

|  |           |
|--|-----------|
| SureSelect Max Library Composition                           | <b>40</b> |
| SureSelect Max UDI Primers Information                       | <b>40</b> |
| Index Primer Pair Strip Tube and Plate Maps                  | <b>41</b> |
| SureSelect Max Index Sequences                               | <b>44</b> |
| Troubleshooting Guide  | <b>52</b> |
| Quick Reference Protocol: Library Prep for Target Enrichment | <b>54</b> |
| Quick Reference Protocol: Library Prep for Whole Genome NGS  | <b>55</b> |

# 1

## Before You Begin

Overview of the Workflow [8](#)  
SureSelect Max Modules Used in the Workflow [9](#)  
Additional Materials Used in the Workflow [10](#)  
Procedural and Safety Notes [12](#)

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

### NOTE

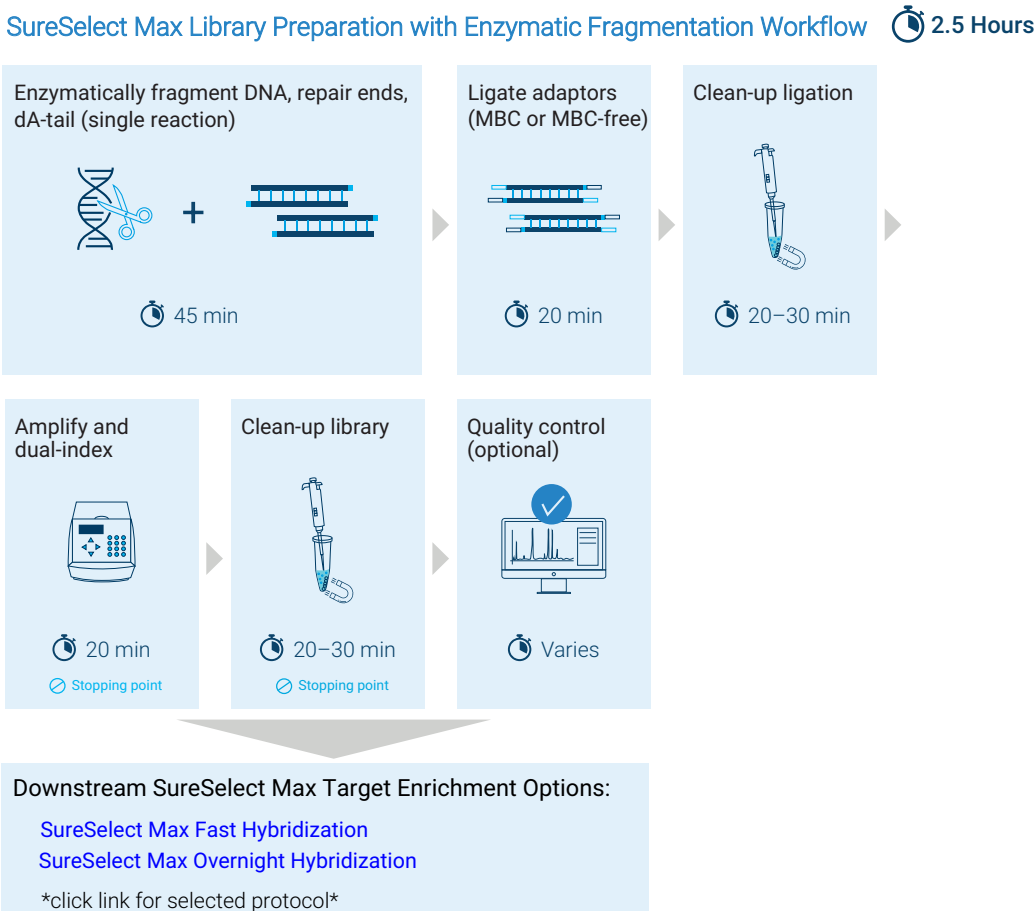
Agilent guarantees performance and provides technical support for the SureSelect Max reagents required for this workflow only when used as directed in this Protocol.

---

# Overview of the Workflow

The SureSelect Max system offers flexible workflow options for preparation and target enrichment of DNA or RNA libraries for NGS, with kits for specific workflow segments in a modular format. This publication provides optimized protocols for the DNA library preparation workflow segment, using enzymatic DNA fragmentation, summarized in [Figure 1](#). The SureSelect Max enzymatic fragmentation module uses a streamlined protocol with DNA fragmentation, end-repair and dA-tailing in a single reaction. For additional flexibility, libraries can be constructed to include or exclude duplex molecular barcodes (MBCs) by using different library adaptors and can be indexed using 384 unique dual indexing (UDI) primers. For detailed protocols see [“DNA Library Preparation with Enzymatic Fragmentation Protocol for Target Enrichment”](#) on page 13. Protocols for downstream target enrichment steps are provided in separate publications.

This publication also supports use of the SureSelect Max Enzymatic Fragmentation Module to prepare DNA libraries for whole genome sequencing using a modified protocol detailed in [“Appendix: Whole Genome Library Preparation Protocol”](#) on page 27.



**Figure 1** Summary of SureSelect Max DNA library preparation using enzymatic fragmentation workflow for downstream target enrichment. The estimated time requirements and optional stopping points are provided in this diagram for reference. Estimates are guidelines for 16 reaction runs using 200 ng high-quality input DNA. Timing for runs using different parameters may vary.

The SureSelect Max system features several improvements over earlier SureSelect platforms:

- Streamlined enzymatic fragmentation library prep protocol with 1-step DNA fragmentation, end-repair and dA-tailing, enhanced chemistry, and support for lower concentration samples
- Enhanced amplification chemistry and master mixed reagents
- Optional pre-capture QC, with support for capture of undiluted library samples
- Enhanced Fast Hyb chemistry and streamlined capture process
- Faster overall turnaround time with shorter, simplified protocol steps

**SureSelect Max reagents are not interchangeable with SureSelect XT HS2 or other SureSelect system reagents.**

## SureSelect Max Modules Used in the Workflow

This publication provides optimized protocols for the library preparation workflow segment, using enzymatic DNA fragmentation. Agilent’s SureSelect reagents required to complete the protocols are summarized in [Table 1](#).

**Table 1** SureSelect Max Kits Used in the DNA Library Prep with Enzymatic Fragmentation Workflow

| Module Description  | 16 Reaction Kits* | 96 Reaction Kits† |
|---|-------------------|-------------------|
| SureSelect Max Enzymatic Fragmentation Library Prep Kit           | G9660A            | G9660B            |
| SureSelect Max Adaptors and UDI Primers Kit for ILM (Select One): |                   |                   |
| MBC Adaptors and UDI Primers 1-16                                 | G9667A            |                   |
| MBC Adaptors and UDI Primers 17-32                                | G9667B            |                   |
| MBC Adaptors and UDI Primers 1-96                                 |                   | G9668A            |
| MBC Adaptors and UDI Primers 97-192                               |                   | G9668B            |
| MBC Adaptors and UDI Primers 193-288                              |                   | G9668C            |
| MBC Adaptors and UDI Primers 289-384                              |                   | G9668D            |
| MBC-Free Adaptors and UDI Primers 1-16                            | G9669A            |                   |
| MBC-Free Adaptors and UDI Primers 17-32                           | G9669B            |                   |
| MBC-Free Adaptors and UDI Primers 1-96                            |                   | G9673A            |
| MBC-Free Adaptors and UDI Primers 97-192                          |                   | G9673B            |
| MBC-Free Adaptors and UDI Primers 193-288                         |                   | G9673C            |
| MBC-Free Adaptors and UDI Primers 289-384                         |                   | G9673D            |
| SureSelect Max Purification Beads                                 | G9962A (5 mL)     | G9962B (30 mL)    |

\* 16-reaction kits contain enough reagents for 2 runs containing 8 samples per run.

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

### NOTE

AMPure XP Beads can replace SureSelect Max Purification Beads in the protocols provided in this user guide (see [Table 2](#) on page 10 for ordering information).

Both types of magnetic purification beads have similar overall performance. The bead types are not, however, identical in composition and the average fragment length and yield results may differ slightly. Use a single bead source in validated protocols.

## Additional Materials Used in the Workflow

See [Table 2](#) through [Table 4](#) for additional reagents and equipment used in the workflow.

**Table 2** Ordering Information for Additional Reagents and Equipment

| Description  | Vendor and Part Number  | Usage Notes   |
|--|---|---|
| gDNA isolation and qualification systems   | Select from <a href="#">Table 3</a> on page 11                            | Select the preparation and qualification systems appropriate for your sample type.  |
| Nucleic acid analysis system   | Select from <a href="#">Table 4</a> on page 11                            | Prepared library QC is optional prior to downstream target enrichment (see <a href="#">page 8</a> ). May also be used for FFPE sample qualification.                                    |
| Qubit BR dsDNA Assay Kit, 100 assays   | Thermo Fisher Scientific p/n Q32850                                       | Use with Thermo Fisher Scientific's Qubit Fluorometer/Assay Tubes (p/n Q33238/Q32856)   |
| 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<br>Tube cap strips | Consult the thermal cycler manufacturer's recommendations                 | —   |
| 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, or equivalent | Requires adapter, p/n C1000-ADAPT, for use with strip tubes   |
| Multichannel and single channel pipettes   | Rainin Pipet-Lite Multi Pipette or equivalent                             | —   |
| Sterile, nuclease-free aerosol barrier pipette tips, vortex mixer, ice bucket, and powder-free gloves            | General laboratory supplier   | —   |
| Magnetic separator   | Thermo Fisher Scientific p/n 12331D or equivalent                         | 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 ring formation.   |
| Nuclease-free Water  | Thermo Fisher Scientific p/n AM9930                                       | Water should not be DEPC-treated.   |
| 100% Ethanol (Ethyl Alcohol, 200 proof)  | Millipore p/n EX0276  | —   |
| Optional: 1X Low TE Buffer (10 mM Tris-HCl, pH 8.0, 0.1 mM EDTA)   | Thermo Fisher Scientific p/n 12090-015, or equivalent                     | Optional solvent for gDNA sample preparation and dilution. Do <b>not</b> use standard TE buffer containing >0.1 mM EDTA; higher EDTA concentrations inhibit the fragmentation reaction. |
| Optional: AMPure XP Kit (5 mL)   | Beckman Coulter Genomics p/n A63880                                       | Optional alternative to SureSelect Max Purification Beads (See <a href="#">Table 1</a> )  |

**Table 3** Recommended DNA Sample Isolation and Qualification Systems

| Description  | Vendor and Part Number | Usage Notes  |
|--|------------------------|--|
| <b>For preparation of high-quality DNA samples</b>               |                        |  |
| QIAamp DNA Mini Kit  | Qiagen                 | Recommended reagents for high-quality DNA sample preparation prior to library preparation.   |
| 50 Samples   | p/n 51304              |  |
| 250 Samples  | p/n 51306              |  |
| <b>For preparation of FFPE DNA samples</b>                       |                        |  |
| QIAamp DNA FFPE Tissue Kit, 50 Samples                           | Qiagen p/n 56404       | Recommended reagents for FFPE gDNA sample preparation prior to library preparation.  |
| Deparaffinization Solution                                       | Qiagen p/n 19093       |  |
| FFPE DNA integrity assessment system:<br>Agilent NGS FFPE QC Kit | Agilent                | Recommended systems for FFPE gDNA qualification prior to library preparation. See <a href="#">Table 4</a> for Agilent TapeStation instrument and accessory ordering information. |
| 16 reactions   | p/n G9700A             |  |
| 96 reactions   | p/n G9700B             |  |
| OR   |                        |  |
| TapeStation Genomic DNA Analysis Consumables:                    | Agilent                |  |
| Genomic DNA ScreenTape   | p/n 5067-5365          |  |
| Genomic DNA Reagents   | p/n 5067-5366          |  |

**Table 4** Recommended Nucleic Acid Analysis Systems

| Analysis System   | Vendor and Part Number Information  | Usage Notes  |
|---|-------------------------------------|--|
| <b>For optional library QC for Target Enrichment workflow</b> |                                     |  |
| Agilent 4200/4150 TapeStation Instrument                      | Agilent p/n G2991AA/G2992AA         | Recommended systems for optional QC of libraries prior to Target Enrichment. (Libraries may also be analyzed using the Agilent 2100 Bioanalyzer instrument and DNA 1000 Kit, p/n 5067-1504.)   |
| Consumables:  |                                     |  |
| 96-well sample plates   | p/n 5042-8502                       |  |
| 96-well plate foil seals                                      | p/n 5067-5154                       |  |
| 8-well tube strips  | p/n 401428                          |  |
| 8-well tube strip caps  | p/n 401425                          |  |
| D1000 ScreenTape  | p/n 5067-5582                       |  |
| D1000 Reagents  | p/n 5067-5583                       |  |
| Agilent 5200/5300/5400 Fragment Analyzer                      | Agilent p/n M5310AA/M5311AA/M5312AA |  |
| Consumables:  |                                     |  |
| NGS Fragment Kit (1-6000 bp)                                  | p/n DNF-473-0500                    |  |
| <b>For required library QC for Whole Genome NGS workflow</b>  |                                     |  |
| Agilent 4200/4150 TapeStation Instrument                      | Agilent p/n G2991AA/G2992AA         | Select one automated electrophoresis system for required whole genome library qualification. May also be used for whole genome library quantification when libraries amplified using $\geq 5$ PCR cycles (see <a href="#">page 29</a> ). |
| Consumables:  |                                     |  |
| 96-well sample plates   | p/n 5042-8502                       |  |
| 96-well plate foil seals                                      | p/n 5067-5154                       |  |
| 8-well tube strips  | p/n 401428                          |  |
| 8-well tube strip caps  | p/n 401425                          |  |
| High Sensitivity D5000 ScreenTape                             | p/n 5067-5592                       |  |
| High Sensitivity D5000 Reagents                               | p/n 5067-5593                       |  |
| Agilent 5200/5300/5400 Fragment Analyzer                      | Agilent p/n M5310AA/M5311AA/M5312AA |  |
| Consumables:  |                                     |  |
| HS NGS Fragment Kit (1-6000 bp)                               | p/n DNF-474-0500                    |  |
| NGS Library Quantification Kit (qPCR-based)                   | Various suppliers                   | Required for quantification of whole genome libraries amplified using $< 5$ PCR cycles (see <a href="#">page 29</a> ).   |

## Procedural and Safety 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 of the following best-practices is recommended to prevent PCR product and nuclease 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 the surfaces that pose the highest risk of contamination daily using a 10% bleach solution, or equivalent.
  - 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.
- Several reagent solutions used in the SureSelect Max protocols are highly viscous. Follow the mixing instructions provided in the protocols.
- Avoid introducing bubbles into reaction mixtures during mixing steps. Before adding sample vials to the thermal cycler for incubation or PCR steps, verify the absence of bubbles at the bottom of the sample wells. If present, spin samples briefly to release the bubbles.
- For each protocol step that requires removal of tube cap strips, reseal the tubes with a fresh strip of 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.
- In general, follow Biosafety Level 1 (BSL1) safety rules.
- Possible stopping points, where samples may be stored at 4°C or –20°C, are marked in the protocol. Do not subject the samples to multiple freeze/thaw cycles.

### CAUTION

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

---

## 2 DNA Library Preparation with Enzymatic Fragmentation Protocol for Target Enrichment

- Step 1. Prepare and qualify genomic DNA samples 15
- Step 2. Fragment, end-repair, and 3'-dA-tail the DNA (Frag/A-Tail) 17
- Step 3. Ligate the adaptor 19
- Step 4. Purify libraries using magnetic purification beads 19
- Step 5. Amplify and index the libraries 21
- Step 6. Purify amplified libraries using magnetic purification beads 22
- Step 7. QC and quantify the libraries (optional) 24

Use the instructions in this section to prepare DNA libraries for downstream target enrichment.

**To prepare DNA libraries for whole genome sequencing, see “Appendix: Whole Genome Library Preparation Protocol” on page 27.**

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. 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.

Library preparation begins with single-reaction enzymatic fragmentation, end-repair and dA-tailing of the DNA fragments. Next, the dA-tailed fragments are ligated to adaptors that either include or exclude duplex molecular barcodes (MBCs). After purification, the library fragments are amplified using unique dual indexing (UDI) primer pairs. After a final purification step, the prepared DNA libraries are ready for later target enrichment and NGS workflow segments. Guidelines are provided at the end of this section for optional QC of the prepared libraries.

For each sample to be sequenced, an individual dual-indexed library is prepared. To process multiple samples together, the protocol includes steps for preparation of reagent mixtures for 8 or 24 samples with overage, followed by distribution to individual DNA samples.

This workflow segment uses the components listed in [Table 5](#). Remove the listed reagents from cold storage and prepare as directed just before use (see the *Where Used* column).

**Table 5** Reagents thawed before use in protocol

| Storage Location   | Kit Component  | Preparative Steps   | Where Used  |
|--|--|---|---|
| SureSelect Max Enzymatic Fragmentation Library Preparation Module, stored at -20°C | Frag/A-Tail Buffer (yellow cap)  | Thaw on ice then keep on ice, vortex to mix.  | <a href="#">page 18</a>                             |
|  | Frag/A-Tail Enzyme Mix (green cap)   | Place on ice just before use. Mix thoroughly by inversion 10X.                          | <a href="#">page 18</a>                             |
|  | Ligation Master Mix (blue cap or bottle)   | Thaw on ice then keep on ice. Mix thoroughly by inversion 10X.                          | <a href="#">page 19</a>                             |
|  | Amplification Master Mix (red cap or bottle)   | Thaw on ice then keep on ice. Mix thoroughly by inversion 10X. Do not vortex.           | <a href="#">page 22</a>                             |
| SureSelect Max Adaptors and UDI Primers Kit for ILM, stored at -20°C               | <b>For MBC-tagged libraries:</b> SureSelect Max MBC Adaptor Oligo Mix, ILM (white cap)<br>OR<br><b>For MBC-free libraries:</b> SureSelect Max MBC-Free Adaptor Oligo Mix (black cap)   | Thaw on ice then keep on ice, vortex to mix.  | <a href="#">page 19</a>                             |
|  | SureSelect Max UDI Primers for ILM (select the specific set of indexes to be used in the run):<br>Index Pairs 1-8 (blue strip)<br>Index Pairs 9-16 (white strip)<br>Index Pairs 17-24 (black strip)<br>Index Pairs 25-32 (red strip)<br>Index Pairs 1-96 (orange plate)<br>Index Pairs 97-192 (blue plate)<br>Index Pairs 193-288 (green plate)<br>Index Pairs 289-384 (red plate) | Thaw on ice then keep on ice, vortex to mix.  | <a href="#">page 22</a>                             |
| +4°C   | SureSelect Max Purification Beads<br>OR<br>AMPure XP Beads   | Equilibrate at room temperature (RT) for at least 30 minutes before use, vortex to mix. | <a href="#">page 19</a> and <a href="#">page 22</a> |

## Step 1. Prepare and qualify genomic DNA samples

Enzymatic DNA fragmentation conditions have been optimized for DNA samples prepared in either 1X Low TE Buffer or nuclease-free water. DNA samples prepared in nuclease-free water may produce slightly smaller insert sizes compared to samples prepared in 1X Low TE Buffer.

**Do not substitute the 1X Low TE Buffer (10 mM Tris-HCl pH 8.0, 0.1 mM EDTA) with standard TE buffer containing >0.1 mM EDTA;** higher EDTA concentrations inhibit the fragmentation reaction.

### 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.

#### 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.
- 3 Place 10–200 ng of each DNA sample in 40  $\mu$ L of nuclease-free water or 1X Low TE Buffer into wells of a thermal cycler-compatible strip tube or PCR plate and hold on ice.

Additional qualification of DNA samples is not required for DNA derived from fresh biological samples. Proceed to ["Step 2. Fragment, end-repair, and 3'-dA-tail the DNA \(Frag/A-Tail\)"](#) on page 17.

### Preparation and qualification of gDNA from FFPE samples

- 1 Prepare gDNA from FFPE tissue sections using a suitable purification system, such as 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.

Do not substitute Buffer ATE with any buffer containing >0.1 mM EDTA; higher EDTA concentrations inhibit the fragmentation reaction.

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

- 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.

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

**Option 1: Qualification using the Agilent Genomic DNA ScreenTape assay DIN score**

The Agilent TapeStation Genomic DNA ScreenTape assay 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 Analyze a 1- $\mu$ L aliquot of each FFPE gDNA sample using the Genomic DNA ScreenTape assay. Follow the instructions provided in the [assay Quick Guide](#).
- b Consult [Table 6](#) for DIN score-based input DNA input guidelines.

**Table 6** 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.

**Option 2: Qualification using the Agilent NGS FFPE QC Kit**

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.

- c Analyze a 1- $\mu$ L aliquot of each FFPE gDNA sample using the Agilent NGS FFPE QC Kit. Follow the instructions provided in the [kit user manual](#).
- d Use the  $\Delta\Delta$ Cq score-based guidelines below (summarized in [Table 7](#)) to determine the appropriate input DNA quantification method for your sample:

For all samples with  $\Delta\Delta$ Cq DNA integrity score  $\leq 1$  (more intact FFPE DNA samples), use the Qubit-based gDNA concentration to determine volume of input DNA needed for the protocol.

For all samples with  $\Delta\Delta$ Cq DNA integrity score  $> 1$  (less intact FFPE DNA samples), 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 7** DNA input guidelines based on  $\Delta\Delta$ Cq DNA integrity score

| $\Delta\Delta$ Cq Score      | DNA Input Guidelines   |
|------------------------------|--|
| $\Delta\Delta$ Cq $\leq 1$ * | 10 ng to 200 ng DNA, based on Qubit Assay quantification         |
| $\Delta\Delta$ Cq $> 1$      | 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.

- 4 After qualification using either method, place 10–200 ng of each FFPE DNA sample in 40 µL of nuclease-free water or 1X Low TE Buffer into wells of a thermal cycler-compatible strip tube or PCR plate and hold on ice.

Proceed to “[Step 2. Fragment, end-repair, and 3'-dA-tail the DNA \(Frag/A-Tail\)](#)” below.

## Step 2. Fragment, end-repair, and 3'-dA-tail the DNA (Frag/A-Tail)

### CAUTION

The SureSelect Max enzymatic fragmentation and library preparation protocols use different reagents and conditions compared to earlier SureSelect protocols. Adhere to the instructions detailed in this publication, including thermal cycler programs and reagent mixing instructions, which differ from those used with earlier SureSelect platforms.

- 1 Preprogram a thermal cycler as shown in [Table 8](#). Set the heated lid to 105°C. Hold at 4°C until samples are added in [step 4](#) on [page 18](#).

**Table 8** Thermal cycler program for Frag/A-Tail reaction (50 µL vol)

| Segment | Temperature | Time   |
|---------|-------------|--|
| 1       | 4°C         | Hold   |
| 2       | 37°C        | Select fragmentation duration based on DNA quality and intended NGS read length (see <a href="#">Table 9</a> ) |
| 3       | 65°C        | 30 minutes   |
| 4       | 4°C         | Hold   |

**Table 9** Fragmentation duration based on DNA quality and NGS read length

| Input type | NGS read length requirement | Fragmentation duration (step 2 of PCR program) |
|------------|-----------------------------|--|
| Intact DNA | 2 ×100 reads                | 20 minutes                                     |
|            | 2 ×150 reads                | 10 minutes                                     |
| FFPE DNA   | 2 ×100 OR 2 ×150 reads      | 15 minutes                                     |

- 2 Prepare the appropriate volume of Frag/A-Tail master mix using the steps below:
  - a Vortex the thawed Frag/A-Tail Buffer at high speed for 5–10 seconds. Spin briefly and keep on ice.
  - b Invert the thawed Frag/A-Tail Enzyme Mix 10 times to homogenize. Spin briefly and keep on ice.
  - c Combine the volume of each reagent listed in [Table 10](#) in a 1.5-mL tube. Seal the tube and mix well by vortexing at high speed for 5 seconds. Spin briefly and keep on ice.

**Table 10** Preparation of Frag/A-Tail master mix

| Reagent                            | Volume for 1 reaction | Volume for 8 reactions (includes excess)* | Volume for 24 reactions (includes excess)† |
|------------------------------------|-----------------------|---|--|
| Frag/A-Tail Buffer (yellow cap)    | 4 µL                  | 36 µL                                     | 104 µL                                     |
| Frag/A-Tail Enzyme Mix (green cap) | 6 µL                  | 54 µL                                     | 156 µL                                     |
| <b>Total</b>                       | <b>10 µL</b>          | <b>90 µL</b>                              | <b>260 µL</b>                              |

\* 16-reaction kits contain enough reagents for 2 runs of 8 samples each using the indicated excess volume. Reagents may be depleted after fewer than 16 reactions when using smaller run sizes.

† 96-reaction kits contain enough reagents for 4 runs of 24 samples each using the indicated excess volume. Reagents may be depleted after fewer than 96 reactions when using smaller run sizes.

- 3 Keep the DNA sample strip or plate on ice while adding 10 µL of the Frag/A-Tail master mix to each sample well containing 40 µL of DNA. Mix by pipetting up and down 15–20 times using a pipette set to 40 µL or cap the wells and vortex at high speed for 5–10 seconds.
- 4 Briefly spin the samples, then immediately place the plate or strip tube in the thermal cycler held at 4°C. Resume the thermal cycling program in [Table 8](#), advancing to the 37°C fragmentation step.

**NOTE**

Remove the magnetic purification beads from cold storage and equilibrate to room temperature (RT) for use on [page 19](#). Keep beads at RT for least 30 minutes before use; beads can be kept at RT through the final pre-capture library purification step on [page 22](#).

## Step 3. Ligate the adaptor

### CAUTION

The Ligation Master Mix used in this step is viscous and requires thorough mixing and careful pipetting. Make sure to follow the mixing instructions below.

Pipette the liquid slowly, ensuring that the full volume is aspirated and dispensed. After addition to the reaction mix, rinse the master mix tip with the sample solution.

- 1 Once the thermal cycling program in [Table 8](#) reaches the final 4°C Hold step, transfer the samples to ice. Preprogram the cycler as show in [Table 11](#) with heated lid off. Hold at 4°C until samples are added in [step 4](#).

**Table 11** Thermal cycler program for ligation (75 µL vol)

| Segment | Temperature | Time       |
|---------|-------------|------------|
| 1       | 4°C         | Hold       |
| 2       | 20°C        | 15 minutes |
| 3       | 4°C         | Hold       |

- 2 Keeping the samples on ice, add 5 µL of the appropriate Adaptor Oligo Mix:
  - For **MBC-tagged** libraries—5 µL of SureSelect Max MBC Adaptor Oligo Mix, ILM (white cap)
  - For **MBC-free** libraries—5 µL of SureSelect Max MBC-Free Adaptor Oligo Mix (black cap)

Mix by pipetting up and down 10 times using a pipette set to 40 µL or cap the wells and vortex at high speed for 5–10 seconds.

- 3 Invert the thawed Ligation Master Mix (blue cap or bottle) 10 times to homogenize. Spin briefly. Add 20 µL of Ligation Master Mix to each sample well. Mix by pipetting up and down slowly at least 15–20 times using a pipette set to 50 µL.

### NOTE

Make sure to add the Adaptor Oligo Mix and the Ligation Master Mix to the samples in separate addition steps, mixing after each addition, as directed above. Adding these components together may increase adaptor-dimer formation and decrease kit performance.

- 4 Briefly spin the samples, then immediately place the plate or strip tube in the thermal cycler and resume the program in [Table 11](#), advancing to the 20°C ligation step.

## Step 4. Purify libraries using magnetic purification beads

### CAUTION

The bead volume used at this step differs in various SureSelect system protocols. Adhere to the instructions provided here; do not use protocols provided for other SureSelect kits.

Once the thermal cycler program in [Table 11](#) reaches the final 4°C hold step, purify the libraries using room-temperature (RT) SureSelect Max Purification Beads or AMPure XP Beads.

Critical purification protocol parameters are summarized for experienced users in [Table 12](#).

**Table 12** Magnetic purification bead cleanup parameters after adaptor ligation

| Parameter   | Value                          |
|---|--------------------------------|
| Volume of RT purification bead suspension added to each sample well | 60 $\mu$ L                     |
| Final elution solvent and volume                                    | 21 $\mu$ L nuclease-free water |
| Amount of eluted sample transferred to fresh well                   | Approximately 20 $\mu$ L       |

- 1 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 all purification steps run on the same day.

- 2 Mix the room-temperature purification beads well until homogeneous and consistent in color.
- 3 Transfer the DNA samples from the thermal cycler to room temperature, then add 60  $\mu$ L of the bead suspension to each sample well.
- 4 Mix by pipetting up and down 15–20 times or cap the wells and vortex at high speed for 5–10 seconds then spin briefly to collect the samples, being careful not to pellet the beads.
- 5 Incubate the bead suspensions 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.
- 8 Continue to keep the plate or strip tube in the magnetic stand while you dispense 200  $\mu$ L of fresh 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](#) and [step 9](#) once for a total of two washes.
- 11 Cap the wells, then briefly spin the samples to collect the residual ethanol. Return the samples to the magnetic stand for 30 seconds. Remove the residual ethanol with a P20 pipette.

**NOTE**

In the drying step below, do not dry the bead pellet to the point that the pellet appears cracked. Elution efficiency is significantly decreased when the bead pellet is excessively dried.

- 12 Dry the samples by keeping the unsealed plate or strip tube at room temperature for 2–5 minutes or until the residual ethanol has just evaporated.
- 13 Elute the library DNA by adding 21  $\mu$ L of nuclease-free water to each sample well.
- 14 Mix by pipetting up and down 10–15 times or cap the wells and vortex at high speed for 5 seconds. Verify that all beads have been resuspended, with no visible clumps in the suspension or bead pellets retained on the sides of the wells. If samples were vortexed, spin briefly to collect the liquid, being careful not to pellet the beads.
- 15 Incubate for 2–5 minutes at room temperature. Use of longer incubation times may improve recovery, especially for longer DNA fragments.
- 16 Put the plate or strip tube in the magnetic stand and leave until the solution is clear (up to 5 minutes).
- 17 Transfer the cleared supernatant (approximately 20  $\mu$ L) to a fresh well. Keep on ice.  
You can discard the beads at this time.

## Step 5. Amplify and index the libraries

- 1 Determine the appropriate SureSelect Max UDI Primer assignment for each sample. See [page 40](#) for information on the UDI primers used to amplify the libraries in this step.

Use a different UDI number for each sample to be sequenced in the same lane.

### NOTE

Agilent's SureSelect 8-bp dual indexes use a uniform numbering system across kit formats and between SureSelect Max for ILM and SureSelect XT HS2 platforms. Do not combine samples indexed with the same-numbered dual index pair from different kit formats for multiplex sequencing.

### CAUTION

The UDI primers are provided in single-use aliquots. To avoid cross-contamination of libraries, do not retain and re-use any residual volume for subsequent experiments.

- 2 Preprogram a thermal cycler as shown in [Table 13](#). Set the heated lid to 105°C. Prewarm the instrument before samples are loaded in [step 6](#) on [page 22](#).

**Table 13** Library indexing/amplification thermal cycler program (50 µL vol)

| Segment | Number of Cycles  | Temperature | Time       |
|---------|---|-------------|------------|
| 1       | 1   | 98°C        | 45 seconds |
| 2       | 7 to 13 based on input DNA quality and quantity (see <a href="#">Table 14</a> ) | 98°C        | 15 seconds |
|         |   | 60°C        | 30 seconds |
|         |   | 72°C        | 30 seconds |
| 3       | 1   | 72°C        | 1 minute   |
| 4       | 1   | 4°C         | Hold       |

**Table 14** Indexing/amplification program cycle number recommendations

| Quality of Input DNA         | Quantity of Input DNA | Cycles    |
|------------------------------|-----------------------|-----------|
| Intact DNA from fresh sample | 200 ng                | 7 cycles  |
|                              | 100 ng                | 8 cycles  |
|                              | 50 ng                 | 9 cycles  |
|                              | 10 ng                 | 10 cycles |
| FFPE sample DNA              | 100 to 200 ng*        | 11 cycles |
|                              | 50 ng*                | 12 cycles |
|                              | 10 ng*                | 13 cycles |

\* For samples qualified by qPCR, use the qPCR-determined quantity of amplifiable DNA to choose cycle number. For samples qualified by DIN value, use the Qubit assay-determined quantity. See [page 15](#) to [page 16](#) for FFPE sample DNA quantification and qualification details.

**NOTE**

To avoid cross-contamination, it is recommended to 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 Mix the thawed Amplification Master Mix (red cap or bottle) thoroughly by 10X inversion then spin briefly.
- 4 Add 25  $\mu\text{L}$  of the Amplification Master Mix to each sample well containing purified DNA library fragments (20  $\mu\text{L}$ ).
- 5 Add 5  $\mu\text{L}$  of the appropriate SureSelect Max UDI primer pair to each reaction.  
Cap 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.
- 6 Place the sample plate or strip tube in the thermal cycler and run the program in [Table 13](#).

**CAUTION**

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

**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.

## Step 6. Purify amplified libraries using magnetic purification beads

Once the thermal cycler program in [Table 13](#) reaches the 4°C hold step, purify the libraries using room-temperature SureSelect Max Purification Beads or AMPure XP Beads.

Critical purification protocol parameters are summarized for experienced users in [Table 15](#).

**Table 15** Magnetic purification bead cleanup parameters after amplification

| Parameter   | Value                                |
|---|--------------------------------------|
| Volume of RT purification bead suspension added to each sample well | 50 $\mu\text{L}$                     |
| Final elution solvent and volume                                    | 15 $\mu\text{L}$ nuclease-free water |
| Amount of eluted sample transferred to fresh well                   | Approximately 14 $\mu\text{L}$       |

- 1 Prepare 400  $\mu\text{L}$  of 70% ethanol per sample, plus excess, for use in [step 8](#).
- 2 Mix the room-temperature purification beads well until homogeneous and consistent in color.
- 3 Transfer the library DNA samples from the thermal cycler to room temperature, then add 50  $\mu\text{L}$  of the bead suspension to each sample well.
- 4 Mix by pipetting up and down 15–20 times or cap the wells and vortex at high speed for 5–10 seconds then spin briefly to collect the samples, being careful not to pellet the beads.
- 5 Incubate the bead suspensions 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 fresh 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](#) and [step 9](#) once for a total of two washes.
- 11 Cap the wells, then briefly spin the samples to collect the residual ethanol. Return the samples to the magnetic stand for 30 seconds. Remove the residual ethanol with a P20 pipette.

**NOTE**

In the drying step below, do not dry the bead pellet to the point that the pellet appears cracked. Elution efficiency is significantly decreased when the bead pellet is excessively dried.

- 12 Dry the samples by keeping the unsealed plate or strip tube at room temperature for 2–5 minutes or until the residual ethanol has just evaporated.
- 13 Elute the library DNA by adding 15  $\mu$ L of nuclease-free water to each sample well.
- 14 Mix by pipetting up and down 10–15 times or cap the wells and vortex at high speed for 5 seconds. Verify that all beads have been resuspended, with no visible clumps in the suspension or bead pellets retained on the sides of the wells. If samples were vortexed, spin briefly to collect the liquid, being careful not to pellet the beads.
- 15 Incubate for 2–5 minutes at room temperature. Use of longer incubation times may improve recovery, especially for longer DNA fragments.
- 16 Put the plate or strip tube in the magnetic stand and leave until the solution is clear (up to 5 minutes).
- 17 Transfer the cleared supernatant (approximately 14  $\mu$ L) to a fresh well. Keep on ice.

You can discard the beads at this time.

**Stopping Point**

If you do not plan to continue to the hybridization workflow segment on same day, seal the wells and store at 4°C overnight or at –20°C for prolonged storage. Remove an aliquot for QC analysis before storage, if appropriate.

## Step 7. QC and quantify the libraries (optional)

QC of the prepared libraries is optional, but quantification is required for hybridization workflows using library normalization, including pre-capture pooling. When normalization is not required, the SureSelect Max Target Enrichment post-capture pooling workflows (both Max Fast Hyb and Max Overnight Hyb) support the use of up to 12  $\mu\text{L}$  of unquantified library samples in hybridization.

For workflows that include pre-capture QC, analyze a sample of each library using one of the platforms in [Table 16](#). Follow the instructions in the linked user guide provided for each assay.

**Table 16** Library analysis options

| Analysis platform                               | Assay used at this step      | Link to assay instructions                                     | Amount of library sample to analyze   |
|---|------------------------------|--|---------------------------------------|
| Agilent 4200/4150 TapeStation system            | D1000 ScreenTape             | <a href="#">Agilent D1000 Assay Quick Guide</a>                | 1 $\mu\text{L}$ of five-fold dilution |
| Agilent 5200/5300/5400 Fragment Analyzer system | NGS Fragment Kit (1-6000 bp) | <a href="#">Agilent NGS Fragment Kit (1-6000 bp) Kit Guide</a> | 2 $\mu\text{L}$ of five-fold dilution |

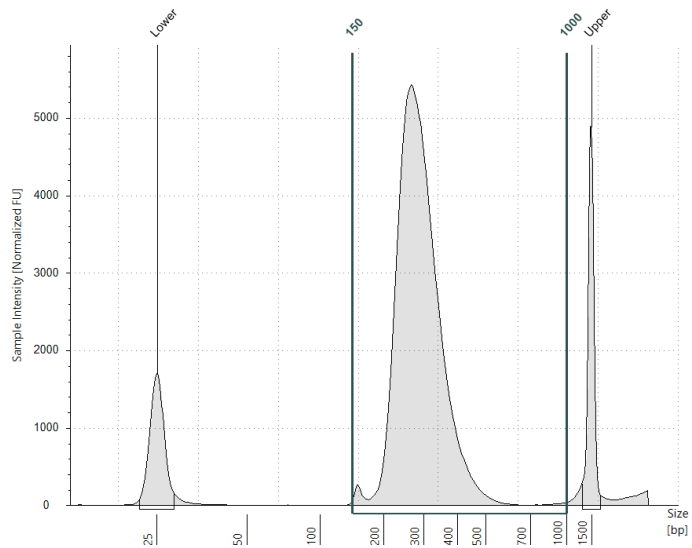
Each analysis method provides an electropherogram showing the size distribution of fragments in the sample and tools for determining the concentration of DNA in the sample. See [Table 17](#) for fragment size distribution guidelines. Representative electropherograms generated using the TapeStation system are provided in [Figure 2](#) and [Figure 3](#) to illustrate typical results.

Using the 150 bp to 1000 bp region of the electropherogram, determine the average fragment size and the library DNA concentration.

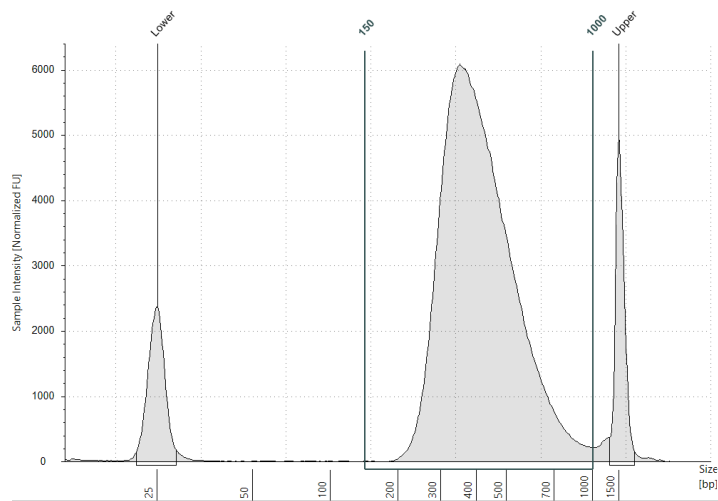
**Table 17** Expected library fragment size guidelines

| Input type | NGS read length used to select fragmentation duration | Expected average fragment size (150–1000 bp region) |
|------------|---|---|
| Intact DNA | 2 $\times$ 100 reads                                  | 250 to 450 bp                                       |
|            | 2 $\times$ 150 reads                                  | 280 to 480 bp                                       |
| FFPE DNA   | 2 $\times$ 100 OR 2 $\times$ 150 reads                | 250 to 360 bp                                       |

Observation of a low molecular weight peak, in addition to the expected library fragment 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, as shown in [Figure 2](#). See *Troubleshooting* on [page 53](#) for additional considerations.



**Figure 2** Library prepared from an FFPE gDNA sample analyzed using a D1000 ScreenTape assay.



**Figure 3** Library prepared from a high-quality gDNA sample analyzed using a D1000 ScreenTape assay.

**Stopping Point** If you do not continue to the hybridization workflow segment on same day, seal the sample wells and store at 4°C overnight or at -20°C for prolonged storage.

The prepared DNA library fragments are ready for target enrichment using the selected workflow option. Proceed to the appropriate SureSelect Max Target Enrichment Module user guide listed in [Table 18](#).

**Table 18** Target Enrichment workflow options

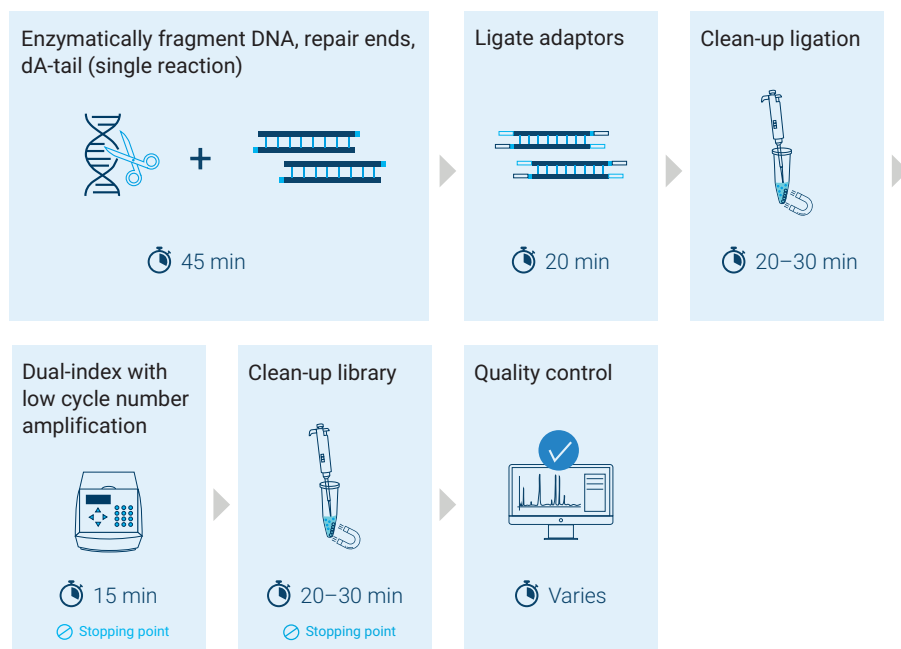
| <b>Workflow option</b>   | <b>Module User Guide link</b> |
|--|-------------------------------|
| Max Fast Hybridization (with pre-capture or post-capture pooling)      | <a href="#">G9689-90000</a>   |
| Max Overnight Hybridization (with pre-capture or post-capture pooling) | <a href="#">G9690-90000</a>   |

### 3 Appendix: Whole Genome Library Preparation Protocol

- Protocol Overview [28](#)
- Step 1. Prepare and qualify genomic DNA samples [28](#)
- Step 2. Fragment, end-repair, and 3'-dA-tail the DNA (Frag/A-Tail) [28](#)
- Step 3. Ligate the adaptor [29](#)
- Step 4. Purify libraries using magnetic purification beads [29](#)
- Step 5. Index and amplify the libraries [29](#)
- Step 6. Purify amplified libraries using magnetic purification beads [30](#)
- Step 7. QC and quantify the libraries [32](#)
- NGS Guidelines [34](#)

This appendix provides an optimized protocol for SureSelect Max Library Preparation for Whole Genome Sequencing using enzymatic DNA fragmentation, summarized in [Figure 4](#). Guidelines are included in this appendix for NGS using the Illumina platform.

**SureSelect Max Library Preparation with Enzymatic Fragmentation Workflow**  **2–2.5 Hours for Whole Genome Sequencing**



**Figure 4** Summary of SureSelect Max DNA library preparation using enzymatic fragmentation for whole genome sequencing. Optional stopping points and estimated time requirements are provided for reference, with timing guidelines based on 16 reaction runs using 200 ng high-quality input DNA. Timing for runs using different protocol parameters may vary.

# Protocol Overview

The whole genome library preparation protocol uses methods and conditions similar to the library preparation protocol for target enrichment provided on [page 14](#) to [page 23](#). This Appendix details the modifications required for whole genome library preparation.

The whole genome library preparation protocol uses the reagents listed in [Table 5](#) on page 14. The SureSelect Max MBC-Free adaptor (black cap) is recommended for use with whole genome sequencing applications. Before you begin, prepare the reagents as directed on [page 14](#).

## Step 1. Prepare and qualify genomic DNA samples

Follow the instructions on [page 15](#) to [page 17](#). Preparation and dilution of gDNA samples using 1X Low TE Buffer is recommended for the whole genome workflow; samples in nuclease-free water can also be used but may produce libraries with smaller average insert sizes.

At the end of this section, you should have samples containing 10–200 ng DNA in 40  $\mu$ L 1X Low TE Buffer (preferred) or nuclease-free water. If you wish to minimize the cycle number in downstream amplification steps, use the maximum input DNA available in the 10–200 ng range.

## Step 2. Fragment, end-repair, and 3'-dA-tail the DNA (Frag/A-Tail)

The Whole Genome Library Preparation protocol uses modified fragmentation conditions to produce longer DNA fragments. After setting up the thermal cycler program below, the instructions on [page 18](#) are applicable to whole genome libraries.

- 1 Preprogram a thermal cycler as shown in [Table 19](#). Set the heated lid to 105°C. Hold at 4°C until samples are loaded.

**Table 19** Thermal cycler program for Frag/A-Tail reaction (50  $\mu$ L vol)

| Step   | Temperature | Time   |
|--------|-------------|--|
| Step 1 | 4°C         | Hold   |
| Step 2 | 37°C        | Select duration based on DNA quality and intended NGS read length ( <a href="#">Table 20</a> ) |
| Step 3 | 65°C        | 30 minutes   |
| Step 4 | 4°C         | Hold   |

**Table 20** Fragmentation duration for whole genome sequencing libraries

| Input type | NGS read length requirement | Fragmentation duration (step 2 of PCR program) |
|------------|-----------------------------|--|
| Intact DNA | 2 $\times$ 150 reads        | 5 minutes                                      |
|            | 2 $\times$ 250 reads        | 2 minutes                                      |
| FFPE DNA   | 2 $\times$ 150 reads        | 5 minutes*                                     |

\* For higher library complexity, fragmentation incubation time can be increased to up to 15 minutes for FFPE samples. Base coverage may be adversely affected when using this modification.

- 2 Proceed to [page 18](#) and follow the instructions in [step 2](#) through [step 4](#) for single-tube fragmentation, end-repair and dA-tailing, using the thermal cycler program in [Table 19](#). At the end of this section, the 50  $\mu$ L DNA samples are held in the thermal cycler.

### Step 3. Ligate the adaptor

Follow the instructions on [page 19](#) ([step 1](#) through [step 4](#)). To prepare MBC-Free whole genome libraries, use the SureSelect Max MBC-Free Adaptor Oligo Mix (black cap) at this step. At the end of this section, the 75  $\mu$ L adaptor-ligated DNA samples are held in the thermal cycler.

### Step 4. Purify libraries using magnetic purification beads

Follow the instructions on [page 20](#) ([step 1](#) through [step 17](#)). At the end of this section, the purified adaptor-ligated DNA samples are in approximately 20  $\mu$ L of nuclease-free water, held on ice.

### Step 5. Index and amplify the libraries

The Whole Genome Library Preparation protocol uses a low PCR cycle number to index and amplify the libraries. Follow the instructions provided below for this step.

- 1 Determine the appropriate SureSelect Max UDI Primer assignment for each sample. Use a different UDI number for each sample to be sequenced in the same lane.
- 2 Preprogram a thermal cycler as shown in [Table 21](#). Set the heated lid to 105°C. Prewarm the instrument before samples are loaded.

#### NOTE

Libraries prepared from 50 ng to 200 ng of intact DNA can be indexed using as few as 3 PCR cycles (see [Table 22](#)). When indexing using <5 PCR cycles, the final libraries must be quantified for sequencing using qPCR. For libraries amplified with  $\geq 5$  cycles, automated electrophoresis may be used for quantification. See [page 32](#) for complete QC guidelines.

**Table 21** Library indexing/amplification thermal cycler program (50  $\mu$ L vol)

| Segment | Number of Cycles   | Temperature | Time       |
|---------|--|-------------|------------|
| 1       | 1  | 98°C        | 45 seconds |
| 2       | 3 to 8 based on input DNA quality and quantity (see <a href="#">Table 22</a> ) | 98°C        | 15 seconds |
|         |  | 60°C        | 30 seconds |
|         |  | 72°C        | 30 seconds |
| 3       | 1  | 72°C        | 1 minute   |
| 4       | 1  | 4°C         | Hold       |

**Table 22** Indexing/amplification program cycle number recommendations

| Quality of Input DNA         | Quantity of Input DNA     | Cycles                  |
|------------------------------|---------------------------|-------------------------|
| Intact DNA from fresh sample | 50 to 200 ng              | 3–5 cycles <sup>*</sup> |
|                              | 10 ng                     | 6 cycles                |
| FFPE sample DNA              | 50 to 200 ng <sup>†</sup> | 7 cycles                |
|                              | 10 ng <sup>†</sup>        | 8 cycles                |

<sup>\*</sup> Libraries amplified using <5 cycles require use of qPCR quantification for sequencing (see [page 32](#)).

<sup>†</sup> For samples qualified by qPCR, use the qPCR-determined quantity of amplifiable DNA to choose cycle number. For samples qualified by DIN value, use the Qubit Assay-determined quantity. See [page 15](#) to [page 16](#) for FFPE sample DNA quantification and qualification details.

- 3 Mix the thawed Amplification Master Mix (red cap or bottle) thoroughly by 10X inversion then spin briefly.
- 4 Add 25 µL of the Amplification Master Mix to each sample well containing purified DNA library fragments (20 µL).
- 5 Add 5 µL of the appropriate SureSelect Max UDI primer pair to each reaction.  
Cap 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.
- 6 Place the sample plate or strip tube in the thermal cycler and run the program in [Table 21](#).

**CAUTION**

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

**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.

## Step 6. Purify amplified libraries using magnetic purification beads

Once the amplification program in [Table 21](#) reaches the final 4°C hold step, move the samples to room temperature and purify using room-temperature SureSelect Max Purification Beads or AMPure XP Beads as directed in this section.

The Whole Genome Library Preparation protocol uses different final purification conditions for 2x250 versus 2x150 read length libraries. These differences and other important purification protocol parameters are summarized in [Table 23](#).

**Table 23** Magnetic purification bead cleanup parameters after amplification

| Step/Parameter  | Value               |  |
|---|---------------------|--|
|   | For 2x150 read NGS  | For 2x250 read NGS                           |
| Two-fold dilution of library prior to purification                  | No                  | Yes (add 50 µL nuclease-free water per well) |
| Volume of RT purification bead suspension added to each sample well | 50 µL               | 60 µL  |
| Duration of initial sample/bead binding incubation                  | 10 minutes          | 10 minutes                                   |
| Final elution solvent and volume                                    | 26 µL Low TE Buffer | 26 µL Low TE Buffer                          |
| Amount of eluted sample transferred to fresh well                   | Approximately 25 µL | Approximately 25 µL                          |

**Important: For libraries to be sequenced using 2 x 250 read length only, dilute each sample two-fold before starting the procedure below.** Dilute by adding 50 µL of nuclease-free water to each 50 µL library sample well.

- 1 Prepare 400 µL of 70% ethanol per sample, plus excess, for use in [step 8](#).
- 2 Mix the room-temperature purification beads well until homogeneous and consistent in color.
- 3 Add the appropriate volume of bead suspension to each sample well:
  - For 2 x 150 NGS libraries: add 50 µL of beads to the 50 µL library samples
  - For 2 x 250 NGS libraries: add 60 µL of beads to the 100 µL diluted library samples
- 4 Mix by pipetting up and down 15–20 times or cap the wells and vortex at high speed for 5–10 seconds then spin briefly to collect the samples, being careful not to pellet the beads.

- 5 Incubate the bead suspensions for 10 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 fresh 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](#) and [step 9](#) once for a total of two washes.
- 11 Cap the wells, then briefly spin the samples to collect the residual ethanol. Return the samples to the magnetic stand for 30 seconds. Remove the residual ethanol with a P20 pipette.

**NOTE**

In the drying step below, do not dry the bead pellet to the point that the pellet appears cracked. Elution efficiency is significantly decreased when the bead pellet is excessively dried.

- 12 Dry the samples by keeping the unsealed plate or strip tube at room temperature for 2–5 minutes or until the residual ethanol has just evaporated.
- 13 Elute the library DNA by adding 26  $\mu$ L of 1X Low TE Buffer to each sample well.
- 14 Mix by pipetting up and down 10–15 times or cap the wells and vortex at high speed for 5 seconds. Verify that all beads have been resuspended, with no visible clumps in the suspension or bead pellets retained on the sides of the wells. If samples were vortexed, spin briefly to collect the liquid, being careful not to pellet the beads.
- 15 Incubate for 2–5 minutes at room temperature. Use of longer incubation times may improve recovery, especially for longer DNA fragments.
- 16 Put the plate or strip tube in the magnetic stand and leave until the solution is clear (up to 5 minutes).
- 17 Transfer the cleared supernatant (approximately 25  $\mu$ L) to a fresh well. Keep on ice. You can discard the beads at this time.

The prepared DNA libraries are ready for NGS after QC and multiplex pooling.

**Stopping Point**

If you do not continue to QC and library pooling on the same day, seal the wells and store at 4°C overnight or at –20°C for prolonged storage.

## Step 7. QC and quantify the libraries

Analyze a sample of each library using the appropriate method(s) summarized in [Table 24](#).

All libraries should be size-qualified using Agilent's TapeStation or Fragment Analyzer system. Libraries indexed using at least 5 PCR cycles (see [Table 22](#) on page 29) can also be quantified using Agilent's automated electrophoresis systems. For libraries prepared using the minimum 3 to 4 amplification cycles, the final libraries must be quantified for sequencing with qPCR.

**Table 24** Library analysis guidelines

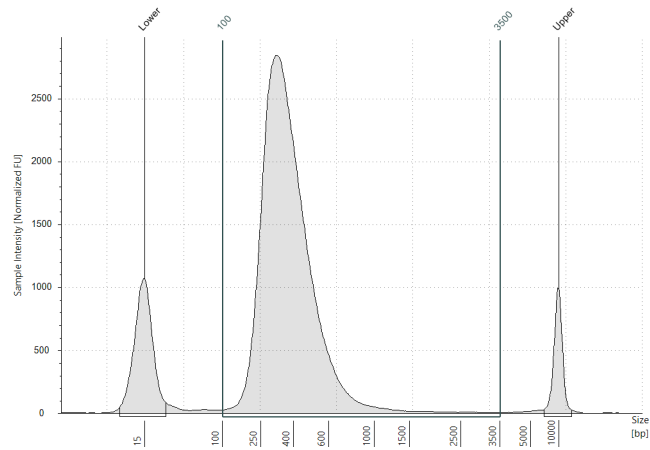
| Purpose  | Analysis platform                               | Assay used at this step                         | Amount of library sample to analyze  |
|--|---|---|--|
| Library fragment size qualification (all libraries) and DNA quantification for libraries amplified using $\geq 5$ cycles | Agilent 4200/4150 TapeStation system            | Agilent High Sensitivity D5000 ScreenTape Assay | 10 ng DNA input libraries: 2 $\mu$ L of 5-fold dilution<br>200 ng DNA input libraries: 2 $\mu$ L of 50-fold dilution |
|  | Agilent 5200/5300/5400 Fragment Analyzer system | Agilent HS NGS Fragment Kit (1-6000 bp)         | 10 ng DNA input libraries: 2 $\mu$ L of 5-fold dilution<br>200 ng DNA input libraries: 2 $\mu$ L of 50-fold dilution |
| DNA quantification for libraries amplified using $<5$ cycles   | qPCR system                                     | qPCR-based NGS Library Quantification Kit       | See manufacturer's protocol  |

Using the 100 bp to 3500 bp region of the electropherogram, determine the average fragment size and the library DNA concentration. See [Table 25](#) for fragment size distribution guidelines. Representative electropherograms generated using the TapeStation system are provided in [Figure 5](#) and [Figure 6](#) to illustrate typical results.

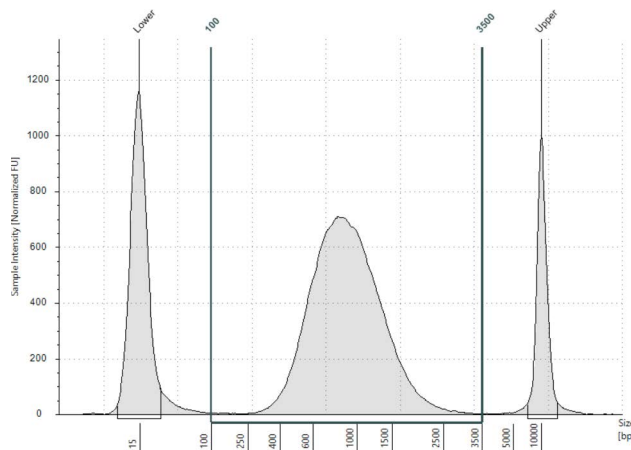
**Table 25** Expected library fragment size guidelines

| Input type | NGS read length used to select fragmentation duration | Expected average fragment size (100–3500 bp region) |
|------------|---|---|
| Intact DNA | 2 $\times$ 150 reads                                  | 510 to 610 bp                                       |
|            | 2 $\times$ 250 reads                                  | 800 to 1000 bp                                      |
| FFPE DNA   | 2 $\times$ 150 reads                                  | 250 to 450 bp                                       |

Quantify libraries prepared with  $<5$  amplification cycles using the appropriate qPCR-based NGS Library Quantification Kit. The approximate library concentration from the automated electrophoresis assay can be used to determine the appropriate dilution scheme for qPCR quantification.



**Figure 5** Library prepared from an FFPE gDNA sample fragmented using 2x150 read length conditions analyzed using a High Sensitivity D5000 ScreenTape assay.



**Figure 6** Library prepared from a high-quality gDNA sample fragmented using 2x250 read length conditions analyzed using a High Sensitivity D5000 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.

# NGS Guidelines

## Library pooling guidelines for multiplex NGS

The SureSelect Max whole genome libraries are ready for pooling for multiplex NGS.

### NOTE

SureSelect Max UDI strip and plate layouts are designed to provide the proper color balance for Illumina's two-channel and four-channel systems. A minimum plexity of four is recommended to ensure that library pools are color balanced. Pools containing any four consecutive SureSelect Max UDIs meet Illumina's guidance for optimal color balance and sequencing performance. Consult Illumina's guidelines for additional color balance and pooling strategy information including two-plex or three-plex pooling considerations.

The number of indexed libraries that may be multiplexed in a single sequencing lane is determined by the output specifications of the sequencer used, together with the amount of sequencing data required per sample 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.

Combine the library samples such that each indexed library is present in equimolar amounts in the NGS pool using one of the following methods. Use the diluent specified by your sequencing provider, such as Low TE, for the dilution steps.

**Method 1:** Dilute each sample to be pooled to the same final concentration (typically 4–15 nM, or the concentration of the most dilute sample) then combine equal volumes of all samples to create the final NGS 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 indexed 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 indexes, and

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

[Table 26](#) shows an example of the amount of 4 indexed samples (of different concentrations) and Low TE needed for a final volume of 20  $\mu\text{L}$  at 10 nM DNA.

**Table 26** Example of volume calculation for total volume of 20  $\mu\text{L}$  at 10 nM concentration

| 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                             |

Store the library pool under the conditions specified by your sequencing provider. Typically, libraries should be stored at  $-20^{\circ}\text{C}$  for short-term storage.

## Sequencing setup and run guidelines

The pooled libraries are ready for 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 (see [page 40](#) for a library structure diagram). Motifs include unique 8-bp P5 and P7 indexes suitable for Illumina sequencing platforms. See [page 40](#) for SureSelect Max UDI information.

Proceed to cluster amplification using the appropriate Illumina Paired-End Cluster Generation Kit. [Table 27](#) provides guidelines for use of several Illumina instrument and chemistry combinations suitable for processing SureSelect Max DNA whole genome NGS libraries quantified using an Agilent automated electrophoresis system. Consult Illumina's documentation for sequencing setup guidelines for qPCR-quantified libraries and for runs using other Illumina instruments.

**Table 27** Illumina Kit Configuration Selection Guidelines

| Instrument           | Run Type                  | Read Length                               | SBS Kit Configuration             | Chemistry    | Seeding Concentration |
|----------------------|---------------------------|---|-----------------------------------|--------------|-----------------------|
| MiSeq                | All Runs                  | 2 $\times$ 150 bp or<br>2 $\times$ 250 bp | 300 Cycle Kit or<br>500 Cycle Kit | v2           | 9–10 pM               |
|                      |                           |   | 600 Cycle Kit                     | v3           | 12–16 pM              |
| NextSeq<br>1000/2000 | All Runs                  | 2 $\times$ 150 bp or<br>2 $\times$ 250 bp | 300 Cycle Kit or<br>600 Cycle Kit | Standard SBS | 650–1000 pM           |
|                      |                           |   | 300 Cycle Kit                     | XLEAP-SBS    | 650–1000 pM           |
| NovaSeq 6000         | Standard<br>Workflow Runs | 2 $\times$ 150 bp or<br>2 $\times$ 250 bp | 300 Cycle Kit or<br>500 Cycle Kit | v1.5         | 300–600 pM            |
|                      | Xp Workflow Runs          | 2 $\times$ 150 bp or<br>2 $\times$ 250 bp | 300 Cycle Kit or<br>500 Cycle Kit | v1.5         | 200–400 pM            |
| iSeq 100             | All Runs                  | 2 $\times$ 150 bp                         | 300 Cycle Kit                     | v2           | 50–150 pM             |
| NextSeq 500/550      | All Runs                  | 2 $\times$ 150 bp                         | 300 Cycle Kit                     | v2.5         | 1.2–1.5 pM            |
| NovaSeq X            | All runs                  | 2 $\times$ 150 bp                         | 300 Cycle Kit                     | v1           | 90–180 pM             |

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 27](#) or provided by Illumina. Follow Illumina's recommendation for a PhiX control in a low-concentration spike-in for improved sequencing quality control.

Set up the sequencing run to generate Read 1 and Read 2 FASTQ files for each sample using the instrument's software in standalone mode or using an Illumina run management tool such as Local Run Manager (LRM), Illumina Experiment Manager (IEM) or BaseSpace. Enter the appropriate **Cycles** or **Read Length** value for your library read length and using 8-bp dual index reads. See [Table 28](#) showing example settings for 2 x 150 bp sequencing.

**Table 28** Run settings for 2 x 150 bp sequencing

| Run Segment  | Cycles/Read Length |
|--------------|--------------------|
| Read 1       | 151*               |
| Index 1 (i7) | 8                  |
| Index 2 (i5) | 8                  |
| Read 2       | 151*               |

\* Follow Illumina's recommendation to add one (1) extra cycle to the desired read length.

Follow Illumina's instructions for each platform and setup software option, incorporating the additional setup guidelines below:

- Each of the sample-level indexes (i7 and i5) requires an 8-bp index read. For complete index sequence information, see [page 44](#) through [page 51](#).
- No custom primers are used for SureSelect Max library sequencing. Leave all *Custom Primers* options for *Read 1*, *Read 2*, *Index 1* and *Index 2* primers cleared/deselected during run setup.
- For runs set up using Illumina's LRM, IEM, or BaseSpace applications, refer to Illumina's instructions and support resources for setting up runs with custom library prep kits and index kits in the selected software. If you need additional SureSelect Max for ILM library sequencing guidance, contact the SureSelect support team (see [page 2](#)) or your local representative.

## Analysis pipeline guidelines

Guidelines are provided below for typical NGS read processing and analysis pipeline steps for SureSelect Max whole genome DNA libraries. Your NGS analysis pipeline may vary.

### MBC-free libraries

- 1 Demultiplexing:** Generate demultiplexed FASTQ files for each sample using Illumina's bcl2fastq, BCL Convert, or a similar software application. This process generates paired-end reads based on the dual indexes and removes sequences with incorrectly paired P5 and P7 indexes.
- 2 Adaptor trimming:** Turn on the adaptor trimming tools in the selected Illumina demultiplexing software to complete adaptor trimming at this step.
- 3 Alignment:** The trimmed reads should be aligned using a suitable tool such as BWA-MEM.

The resulting BAM files are ready for downstream analysis including variant discovery.

## MBC-tagged libraries

- 1 Demultiplexing:** Generate demultiplexed FASTQ files for each sample using Illumina's bcl2fastq, BCL Convert, or a similar software application. This process generates paired-end reads based on the dual indexes and removes sequences with incorrectly paired P5 and P7 indexes.

Turn off the MBC/UMI trimming options in Illumina's demultiplexing software to allow proper adaptor processing and use of the MBCs by downstream tools.

### NOTE

If your libraries are MBC-tagged but your sequence analysis pipeline excludes MBCs, you can remove the first 5 bases from Read 1 and Read 2 by masking or trimming before proceeding to downstream analysis.

If demultiplexing using bcl2fastq, MBCs may be masked by including the base mask **N5Y\*,I8,I8,N5Y\*** (where \* is replaced with the remaining read length after subtracting the 5 masked bases, e.g., use **N5Y146,I8,I8,N5Y146** for 2x150 NGS set up as shown in [Table 28](#) on page 36). The sum of the values following N and Y must match the read length value in the RunInfo.xml file.

If demultiplexing using BCL Convert, MBCs may be trimmed by including the following string in the sample sheet header: **OverrideCycles,N5Y\*;I8,I8,N5Y\*** (where \* is replaced with read length after trimming, e.g., use **N5Y146;I8;I8;N5Y146** for 2x150 NGS set up as shown in [Table 28](#) on page 36). The sum of the values following N and Y must match the read length value in the RunInfo.xml file.

Alternatively, the first 5 bases may be trimmed from the demultiplexed FASTQ files using a suitable processing tool (e.g., fgbio).

- 2 MBC-adaptor processing:** Use a suitable processing tool of your choice to trim and collect inline MBCs from each sequencing read. For example, MBC processing could be conducted with the fgbio best practice consensus pipeline, as described in: <https://github.com/fulcrumgenomics/fgbio/blob/main/docs/best-practice-consensus-pipeline.md>. The output includes trimmed, deduplicated reads in unaligned BAM format with single-stranded MBC consensus reads. For generation of double-stranded MBC consensus reads, refer to <https://fulcrumgenomics.github.io/fgbio/tools/latest/CallDuplexConsensusReads.html>.
  - Inline MBCs are added to both ends of the DNA inserts in the assay. To collect the MBCs and process them for analysis, one approach is to trim off 5 bases at the beginning of each read, take the first 3 of the 5 bases as MBC, and discard the remaining 2.
- 3 Alignment:**
  - DNA library alignment can be completed as part of the fgbio best practice consensus pipeline described at the link above or aligned using another suitable alignment tool of your choice.

The resulting aligned BAM files are ready for downstream analysis including variant discovery.

## 4 Reference

|  |    |
|--|----|
| Reagent Kit Contents   | 39 |
| SureSelect Max Library Composition                           | 40 |
| SureSelect Max UDI Primers Information                       | 40 |
| Troubleshooting Guide  | 52 |
| Quick Reference Protocol: Library Prep for Target Enrichment | 54 |
| Quick Reference Protocol: Library Prep for Whole Genome NGS  | 55 |

This section contains reference information, including Reagent Kit contents, index sequences, troubleshooting information and quick-reference protocols for experienced users.

## Reagent Kit Contents

SureSelect Max DNA Library Preparation with enzymatic fragmentation uses the kits listed in [Table 29](#). Detailed contents of the multi-part component kits are shown in [Table 30](#) through [Table 32](#).

**Table 29** Kits for SureSelect Max DNA Library Preparation with Enzymatic Fragmentation

| Purchased Kit  | Included Component Kits   | Component Kit Part Number                         |  | Storage Condition |
|--|---|---|--|-------------------|
|  |   | 16 Reactions                                      | 96 Reactions   |                   |
| SureSelect Max Enzymatic Fragmentation Library Preparation Kit | SureSelect Max Enzymatic Fragmentation Library Preparation Module | 5280-0063   | 5280-0064  | -20°C             |
| SureSelect Max Adaptors and UDI Primers Kit for ILM            | SureSelect Max MBC Adaptor Oligo Mix for ILM                      | 5282-0124   | 5282-0125  | -20°C             |
|  | <b>OR</b><br>SureSelect Max MBC-Free Adaptor Oligo Mix for ILM    | <b>OR</b><br>5282-0126                            | <b>OR</b><br>5282-0127   |                   |
|  | SureSelect Max UDI Primers for ILM                                | 5282-0138 (Index 1-16)<br>5282-0119 (Index 17-32) | 5282-0120 (Index 1-96)<br>5282-0121 (Index 97-192)<br>5282-0122 (Index 193-288)<br>5282-0123 (Index 289-384) | -20°C             |
| SureSelect Max Purification Beads                              |   | 5282-0225   | 5282-0226  | +4°C              |

## Component Kit Details

**Table 30** SureSelect Max Library Preparation Module content

| Kit Component            | 16 Reaction Kit (p/n 5280-0063) | 96 Reaction Kit (p/n 5280-0064) |
|--------------------------|---------------------------------|---------------------------------|
| Frag/A-Tail Enzyme Mix   | tube with green cap             | tube with green cap             |
| Frag/A-Tail Buffer       | tube with yellow cap            | tube with yellow cap            |
| Ligation Master Mix      | tube with blue cap              | bottle                          |
| Amplification Master Mix | tube with red cap               | bottle                          |

**Table 31** SureSelect Max Adaptor Oligo Mix for ILM options

| Kit Component                                 | 16 Reaction Kits    | 96 Reaction Kits    |
|---|---------------------|---------------------|
| SureSelect MBC Adaptor Oligo Mix for ILM      | tube with white cap | tube with white cap |
| SureSelect MBC-Free Adaptor Oligo Mix for ILM | tube with black cap | tube with black cap |

**Table 32** SureSelect Max UDI Primers for ILM options

| Kit Component                       | 16 Reaction Kit Format   | 96 Reaction Kit Format   |
|-------------------------------------|--|--|
| SureSelect Max UDI Primers for ILM* | Blue 8-well strip tube (index pairs 1-8), AND<br>White 8-well strip tube (index pairs 9-16)<br><b>OR</b><br>Black 8-well strip tube (index pairs 17-24) AND<br>red 8-well strip tube (index pairs 25-32) | Orange 96-well plate (index pairs 1-96), OR<br>Blue 96-well plate (index pairs 97-192), OR<br>Green 96-well plate (index pairs 193-288), OR<br>Red 96-well plate (index pairs 289-384) |

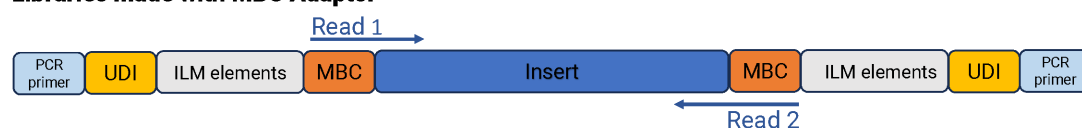
\* See [page 41](#) through [page 43](#) for index strip and plate position maps; see [page 44](#) through [page 51](#) for index pair sequence information.

## SureSelect Max Library Composition

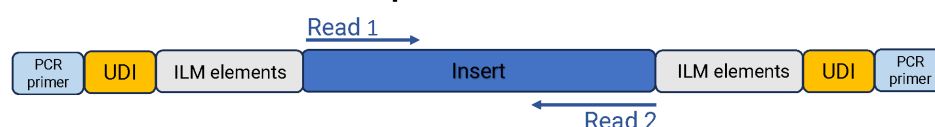
The final SureSelect Max library fragments prepared using Agilent's Illumina-compatible kit modules are shown in [Figure 7](#). Each fragment in the prepared library contains one target insert surrounded by sequence motifs required for multiplexed sequencing using Illumina sequencers. The libraries are ready for sequencing using standard Illumina paired-end primers and chemistry.

Each library DNA fragment contains a unique 8-bp P5 and P7 index suitable for Illumina sequencing platforms. See "[SureSelect Max UDI Primers Information](#)" below for additional information.

### Libraries made with MBC Adaptor



### Libraries made with MBC-Free Adaptor



**Figure 7** Content of SureSelect Max for ILM sequencing libraries. Each fragment contains one target insert (dark blue) surrounded by the following elements: Illumina paired-end sequencing elements (gray), unique dual indexes (gold), library PCR primers (light blue) and optional molecular barcode (MBC) sequences. MBCs include a 3-bp degenerate barcode plus dark bases.

## SureSelect Max UDI Primers Information

The SureSelect Max unique dual indexing (UDI) Primers are provided in pre-combined pairs of indexes. Each member of the primer pair contains a unique 8-bp P5 or P7 index, resulting in dual-indexed NGS libraries. One primer pair is provided in each well of 8-well strip tubes (16 reaction kits; see [Figure 8](#)) or of 96-well plates (96 reaction kits; see [page 42](#) to [page 43](#) for plate maps). Each well contains a single-use aliquot of a specific pair of forward plus reverse primers.

The nucleotide sequence of the index portion of each primer is provided in [Table 37](#) on page 44 through [Table 44](#) on page 51. Index sequences can also be obtained by downloading the [SureSelect Max Index Sequence Resource](#) Excel spreadsheet from Agilent.com.

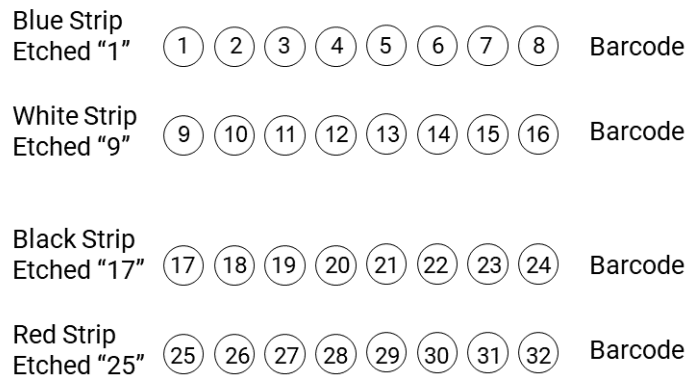
### NOTE

Clicking the Excel spreadsheet link from Agilent.com automatically downloads the index spreadsheet to the default folder for downloaded files saved by your web browser. Locate the file in the folder and open it in Microsoft Excel or another compatible spreadsheet program. The first tab of the spreadsheet provides instructions for use of the spreadsheet contents.

In [Table 37](#) through [Table 44](#) and in the downloadable Excel spreadsheet, P7 indexes are shown in forward orientation, applicable to any of the supported Illumina platforms. P5 indexes are shown in both forward and reverse complement orientations. Refer to Illumina support documentation and resources to determine the correct P5 index orientation for your application.

## Index Primer Pair Strip Tube and Plate Maps

SureSelect Max UDI Primers 1-16 and 17-32 (provided with 16 reaction kits) are supplied in sets of two 8-well strip tubes as detailed below.



**Figure 8** Map of the SureSelect Max UDI Primers for ILM strip tubes in 16 reaction kits

The blue strip contains Index Primer Pairs 1-8, with pair #1 supplied in the well proximal to the numeral **1** etched on the strip's plastic end tab.

The white strip contains Index Primer Pairs 9-16, with pair #9 supplied in the well proximal to the numeral **9** etched on the strip's plastic end tab.

The black strip contains Index Primer Pairs 17-24, with pair #17 supplied in the well proximal to the numeral **17** etched on the strip's plastic end tab.

The red strip contains Index Primer Pairs 25-32, with pair #25 supplied in the well proximal to the numeral **25** etched on the strip's plastic end tab.

When using the strip tube-supplied index primer pairs in the library preparation protocol, pierce the foil seal of the appropriate well with a pipette tip just before pipetting the solution. If the foil seal for any unused wells is disrupted during use, the unused wells may be re-sealed using the provided fresh foil seal strips. The provided foil strips may also be used to re-seal used wells to prevent index pair cross-contamination during subsequent use.

Plate positions of the SureSelect Max UDI Primers for ILM provided with 96 reaction kits are shown in [Table 33](#) through [Table 36](#).

### CAUTION

The SureSelect Max UDI Primers are provided in single-use aliquots. To avoid cross-contamination of libraries, use each well in only one library preparation reaction. Do not retain and re-use any residual volume for subsequent experiments.

**Table 33** Plate map for SureSelect Max UDI Primers 1-96, provided in orange 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 |

**Table 34** Plate map for SureSelect Max UDI Primers 97-192, provided in blue plate

|   | 1   | 2   | 3   | 4   | 5   | 6   | 7   | 8   | 9   | 10  | 11  | 12  |
|---|-----|-----|-----|-----|-----|-----|-----|-----|-----|-----|-----|-----|
| A | 97  | 105 | 113 | 121 | 129 | 137 | 145 | 153 | 161 | 169 | 177 | 185 |
| B | 98  | 106 | 114 | 122 | 130 | 138 | 146 | 154 | 162 | 170 | 178 | 186 |
| C | 99  | 107 | 115 | 123 | 131 | 139 | 147 | 155 | 163 | 171 | 179 | 187 |
| D | 100 | 108 | 116 | 124 | 132 | 140 | 148 | 156 | 164 | 172 | 180 | 188 |
| E | 101 | 109 | 117 | 125 | 133 | 141 | 149 | 157 | 165 | 173 | 181 | 189 |
| F | 102 | 110 | 118 | 126 | 134 | 142 | 150 | 158 | 166 | 174 | 182 | 190 |
| G | 103 | 111 | 119 | 127 | 135 | 143 | 151 | 159 | 167 | 175 | 183 | 191 |
| H | 104 | 112 | 120 | 128 | 136 | 144 | 152 | 160 | 168 | 176 | 184 | 192 |

**Table 35** Plate map for SureSelect Max UDI Primers 193-288, provided in green plate

|   | 1   | 2   | 3   | 4   | 5   | 6   | 7   | 8   | 9   | 10  | 11  | 12  |
|---|-----|-----|-----|-----|-----|-----|-----|-----|-----|-----|-----|-----|
| A | 193 | 201 | 209 | 217 | 225 | 233 | 241 | 249 | 257 | 265 | 273 | 281 |
| B | 194 | 202 | 210 | 218 | 226 | 234 | 242 | 250 | 258 | 266 | 274 | 282 |
| C | 195 | 203 | 211 | 219 | 227 | 235 | 243 | 251 | 259 | 267 | 275 | 283 |
| D | 196 | 204 | 212 | 220 | 228 | 236 | 244 | 252 | 260 | 268 | 276 | 284 |
| E | 197 | 205 | 213 | 221 | 229 | 237 | 245 | 253 | 261 | 269 | 277 | 285 |
| F | 198 | 206 | 214 | 222 | 230 | 238 | 246 | 254 | 262 | 270 | 278 | 286 |
| G | 199 | 207 | 215 | 223 | 231 | 239 | 247 | 255 | 263 | 271 | 279 | 287 |
| H | 200 | 208 | 216 | 224 | 232 | 240 | 248 | 256 | 264 | 272 | 280 | 288 |

**Table 36** Plate map for SureSelect Max UDI Primers 289-384, provided in red plate

|   | 1   | 2   | 3   | 4   | 5   | 6   | 7   | 8   | 9   | 10  | 11  | 12  |
|---|-----|-----|-----|-----|-----|-----|-----|-----|-----|-----|-----|-----|
| A | 289 | 297 | 305 | 313 | 321 | 329 | 337 | 345 | 353 | 361 | 369 | 377 |
| B | 290 | 298 | 306 | 314 | 322 | 330 | 338 | 346 | 354 | 362 | 370 | 378 |
| C | 291 | 299 | 307 | 315 | 323 | 331 | 339 | 347 | 355 | 363 | 371 | 379 |
| D | 292 | 300 | 308 | 316 | 324 | 332 | 340 | 348 | 356 | 364 | 372 | 380 |
| E | 293 | 301 | 309 | 317 | 325 | 333 | 341 | 349 | 357 | 365 | 373 | 381 |
| F | 294 | 302 | 310 | 318 | 326 | 334 | 342 | 350 | 358 | 366 | 374 | 382 |
| G | 295 | 303 | 311 | 319 | 327 | 335 | 343 | 351 | 359 | 367 | 375 | 383 |
| H | 296 | 304 | 312 | 320 | 328 | 336 | 344 | 352 | 360 | 368 | 376 | 384 |

## SureSelect Max Index Sequences

Table 37 SureSelect Max Index Primer Pairs 1–48, provided in orange 96-well plate or in strip tubes

| Primer Pair # | Well | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement | Primer Pair # | Well | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement |
|---------------|------|------------------|------------------|-----------------------------|---------------|------|------------------|------------------|-----------------------------|
| 1             | A01  | CAAGGTGA         | ATGGTTAG         | CTAACCAT                    | 25            | A04  | AGATGGAT         | TGGCACCA         | TGGTGCCA                    |
| 2             | B01  | TAGACCAA         | CAAGGTGA         | TCACCTTG                    | 26            | B04  | GAATTGTG         | AGATGGAT         | ATCCATCT                    |
| 3             | C01  | AGTCGCGA         | TAGACCAA         | TTGGTCTA                    | 27            | C04  | GAGCACTG         | GAATTGTG         | CACAATTC                    |
| 4             | D01  | CGGTAGAG         | AGTCGCGA         | TCGCGACT                    | 28            | D04  | GTTGCGGA         | GAGCACTG         | CAGTGCTC                    |
| 5             | E01  | TCAGCATC         | AAGGAGCG         | CGCTCCTT                    | 29            | E04  | AATGGAAC         | GTTGCGGA         | TCCGCAAC                    |
| 6             | F01  | AGAAGCAA         | TCAGCATC         | GATGCTGA                    | 30            | F04  | TCAGAGGT         | AATGGAAC         | GTTCCATT                    |
| 7             | G01  | GCAGGTTC         | AGAAGCAA         | TTGCTTCT                    | 31            | G04  | GCAACAAT         | TCAGAGGT         | ACCTCTGA                    |
| 8             | H01  | AAGTGTCT         | GCAGGTTC         | GAACCTGC                    | 32            | H04  | GTCGATCG         | GCAACAAT         | ATTGTTGC                    |
| 9             | A02  | CTACCGAA         | AAGTGTCT         | AGACACTT                    | 33            | A05  | ATGGTAGC         | GTCGATCG         | CGATCGAC                    |
| 10            | B02  | TAGAGCTC         | CTACCGAA         | TTCGGTAG                    | 34            | B05  | CGCCAATT         | ATGGTAGC         | GCTACCAT                    |
| 11            | C02  | ATGTCAAG         | TAGAGCTC         | GAGCTCTA                    | 35            | C05  | GACAATTG         | CGCCAATT         | AATTGGCG                    |
| 12            | D02  | GCATCATA         | ATGTCAAG         | CTTGACAT                    | 36            | D05  | ATATTCCG         | GACAATTG         | CAATTGTC                    |
| 13            | E02  | GACTTGAC         | GCATCATA         | TATGATGC                    | 37            | E05  | TCTACCTC         | ATATTCCG         | CGGAATAT                    |
| 14            | F02  | CTACAATG         | GACTTGAC         | GTCAAGTC                    | 38            | F05  | TCGTCGTG         | TCTACCTC         | GAGGTAGA                    |
| 15            | G02  | TCTCAGCA         | CTACAATG         | CATTGTAG                    | 39            | G05  | ATGAGAAC         | TCGTCGTG         | CACGACGA                    |
| 16            | H02  | AGACACAC         | TCTCAGCA         | TGCTGAGA                    | 40            | H05  | GTCCTATA         | ATGAGAAC         | GTTTCAT                     |
| 17            | A03  | CAGGTCTG         | AGACACAC         | GTGTGTCT                    | 41            | A06  | AATGACCA         | GTCCTATA         | TATAGGAC                    |
| 18            | B03  | AATACGCG         | CAGGTCTG         | CAGACCTG                    | 42            | B06  | CAGACGCT         | AATGACCA         | TGGTCATT                    |
| 19            | C03  | GCACACAT         | AATACGCG         | CGCGTATT                    | 43            | C06  | TCGAACTG         | CAGACGCT         | AGCGTCTG                    |
| 20            | D03  | CTTGCATA         | GCACACAT         | ATGTGTGC                    | 44            | D06  | CGCTTCCA         | TCGAACTG         | CAGTTCGA                    |
| 21            | E03  | ATCCTCTT         | CTTGCATA         | TATGCAAG                    | 45            | E06  | TATTCTG          | CGCTTCCA         | TGGAAGCG                    |
| 22            | F03  | GCACCTAA         | ATCCTCTT         | AAGAGGAT                    | 46            | F06  | CAAGTTAC         | TATTCTG          | CAGGAATA                    |
| 23            | G03  | TGCTGCTC         | GCACCTAA         | TTAGGTGC                    | 47            | G06  | CAGAGCAG         | CAAGTTAC         | GTAACCTG                    |
| 24            | H03  | TGGCACCA         | TGCTGCTC         | GAGCAGCA                    | 48            | H06  | CGCGCAAT         | CAGAGCAG         | CTGCTCTG                    |

**Table 38** SureSelect Max Index Primer Pairs 49–96, provided in orange 96-well plate

| Primer Pair # | Well | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement | Primer Pair # | Well | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement |
|---------------|------|------------------|------------------|-----------------------------|---------------|------|------------------|------------------|-----------------------------|
| 49            | A07  | TGAGGAGT         | CGCGCAAT         | ATTGCGCG                    | 73            | A10  | AACGCATT         | ATAGTGAC         | GTCACTAT                    |
| 50            | B07  | ATGACGAA         | TGAGGAGT         | ACTCCTCA                    | 74            | B10  | CAGTTGCG         | AACGCATT         | AATGCGTT                    |
| 51            | C07  | TACGGCGA         | ATGACGAA         | TTCGTCAT                    | 75            | C10  | TGCCTCGA         | CAGTTGCG         | CGCAACTG                    |
| 52            | D07  | AGCGAGTT         | TACGGCGA         | TCGCCGTA                    | 76            | D10  | AAGGCTTA         | TGCCTCGA         | TCGAGGCA                    |
| 53            | E07  | TGTATCAC         | AGCGAGTT         | AACTCGCT                    | 77            | E10  | GCAATGAA         | AAGGCTTA         | TAAGCCTT                    |
| 54            | F07  | GATCGCCT         | TGTATCAC         | GTGATACA                    | 78            | F10  | AAGAACCT         | GCAATGAA         | TTCATTGC                    |
| 55            | G07  | GACTCAAT         | GATCGCCT         | AGGCGATC                    | 79            | G10  | CTGTGCCT         | AAGAACCT         | AGGTTCTT                    |
| 56            | H07  | CAGCTTGC         | GACTCAAT         | ATTGAGTC                    | 80            | H10  | TACGTAGC         | CTGTGCCT         | AGGCACAG                    |
| 57            | A08  | AGCTGAAG         | CAGCTTGC         | GCAAGCTG                    | 81            | A11  | AAGTGGAC         | TACGTAGC         | GCTACGTA                    |
| 58            | B08  | ATTCCGTG         | AGCTGAAG         | CTTCAGCT                    | 82            | B11  | CAACCGTG         | AAGTGGAC         | GTCCACTT                    |
| 59            | C08  | TATGCCGC         | ATTCCGTG         | CACGGAAT                    | 83            | C11  | CTGTTGTT         | CAACCGTG         | CACGGTTG                    |
| 60            | D08  | TCAGCTCA         | TATGCCGC         | GCGGCATA                    | 84            | D11  | GCACGATG         | CTGTTGTT         | AACAACAG                    |
| 61            | E08  | AACTGCAA         | TCAGCTCA         | TGAGCTGA                    | 85            | E11  | GTACGGAC         | GCACGATG         | CATCGTGC                    |
| 62            | F08  | ATTAGGAG         | AACTGCAA         | TTGCAGTT                    | 86            | F11  | CTCCAAGC         | GTACGGAC         | GTCCGTAC                    |
| 63            | G08  | CAGCAATA         | ATTAGGAG         | CTCCTAAT                    | 87            | G11  | TAGTCTGA         | CTCCAAGC         | GCTTGGAG                    |
| 64            | H08  | GCCAAGCT         | CAGCAATA         | TATTGCTG                    | 88            | H11  | TTCGCCGT         | TAGTCTGA         | TCAGACTA                    |
| 65            | A09  | TCCGTTAA         | GCCAAGCT         | AGCTTGGC                    | 89            | A12  | GAACTAAG         | ATACGAAG         | CTTCGTAT                    |
| 66            | B09  | GTGCAACG         | TCCGTTAA         | TTAACGGA                    | 90            | B12  | AAGCCATC         | GAGATTCA         | TGAATCTC                    |
| 67            | C09  | AGTAACGC         | GTGCAACG         | CGTTGCAC                    | 91            | C12  | AACTCTTG         | AAGCCATC         | GATGGCTT                    |
| 68            | D09  | CATAGCCA         | AGTAACGC         | GCGTFACT                    | 92            | D12  | GTAGTCAT         | AACTCTTG         | CAAGAGTT                    |
| 69            | E09  | CACTAGTA         | CATAGCCA         | TGGCTATG                    | 93            | E12  | CTCGCTAG         | GTAGTCAT         | ATGACTAC                    |
| 70            | F09  | TTAGTGCG         | CACTAGTA         | TACTAGTG                    | 94            | F12  | AGTCTTCA         | CAGTATCA         | TGATACTG                    |
| 71            | G09  | TCGATACA         | TTAGTGCG         | CGCACTAA                    | 95            | G12  | TCAAGCTA         | CTTCGTAC         | GTACGAAG                    |
| 72            | H09  | ATAGTGAC         | TCGATACA         | TGTATCGA                    | 96            | H12  | CTTATCCT         | TCAAGCTA         | TAGCTTGA                    |

**Table 39** SureSelect Max Index Primer Pairs 97–144, provided in blue 96-well plate

| Primer Pair # | Well | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement | Primer Pair # | Well | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement |
|---------------|------|------------------|------------------|-----------------------------|---------------|------|------------------|------------------|-----------------------------|
| 97            | A01  | TCATCCTT         | CTTATCCT         | AGGATAAG                    | 121           | A04  | CAGGCAGA         | AGACGCCT         | AGGCGTCT                    |
| 98            | B01  | AACACTCT         | TCATCCTT         | AAGGATGA                    | 122           | B04  | TCCGCGAT         | CAGGCAGA         | TCTGCCTG                    |
| 99            | C01  | CACCTAGA         | AACACTCT         | AGAGTGTT                    | 123           | C04  | CTCGTACG         | TCCGCGAT         | ATCGCGGA                    |
| 100           | D01  | AGTTCATG         | CACCTAGA         | TCTAGGTG                    | 124           | D04  | CACACATA         | CTCGTACG         | CGTACGAG                    |
| 101           | E01  | GTTGGTGT         | AGTTCATG         | CATGAACT                    | 125           | E04  | CGTCAAGA         | CACACATA         | TATGTGTG                    |
| 102           | F01  | GCTACGCA         | GTTGGTGT         | ACACCAAC                    | 126           | F04  | TTCGCGCA         | CGTCAAGA         | TCTTGACG                    |
| 103           | G01  | TCAACTGC         | GCTACGCA         | TGCGTAGC                    | 127           | G04  | CGACTACG         | TTCGCGCA         | TGCGCGAA                    |
| 104           | H01  | AAGCGAAT         | TCAACTGC         | GCAGTTGA                    | 128           | H04  | GAAGGTAT         | CGACTACG         | CGTAGTCG                    |
| 105           | A02  | GTGTTACA         | AAGCGAAT         | ATTCGCTT                    | 129           | A05  | TTGGCATG         | GAAGGTAT         | ATACCTTC                    |
| 106           | B02  | CAAGCCAT         | GTGTTACA         | TGTAACAC                    | 130           | B05  | CGAATTCA         | TTGGCATG         | CATGCCAA                    |
| 107           | C02  | CTCTCGTG         | CAAGCCAT         | ATGGCTTG                    | 131           | C05  | TTAGTTGC         | CGAATTCA         | TGAATTCG                    |
| 108           | D02  | TCGACAAC         | CTCTCGTG         | CACGAGAG                    | 132           | D05  | GATGCCAA         | TTAGTTGC         | GCAACTAA                    |
| 109           | E02  | TCGATGTT         | TCGACAAC         | GTTGTCGA                    | 133           | E05  | AGTTGCCG         | GATGCCAA         | TTGGCATC                    |
| 110           | F02  | CAAGGAAG         | TCGATGTT         | AACATCGA                    | 134           | F05  | GTCCACCT         | AGTTGCCG         | CGGCAACT                    |
| 111           | G02  | ATTGATGC         | AGAGAATC         | GATTCTCT                    | 135           | G05  | ATCAAGGT         | GTCCACCT         | AGGTGGAC                    |
| 112           | H02  | TCGCAGAT         | TTGATGGC         | GCCATCAA                    | 136           | H05  | GAACCAGA         | ATCAAGGT         | ACCTTGAT                    |
| 113           | A03  | GCAGAGAC         | TCGCAGAT         | ATCTGCGA                    | 137           | A06  | CATGTTCT         | GAACCAGA         | TCTGGTTC                    |
| 114           | B03  | CTGCGAGA         | GCAGAGAC         | GTCTCTGC                    | 138           | B06  | TCACTGTG         | CATGTTCT         | AGAACATG                    |
| 115           | C03  | CAACCAAC         | CTGCGAGA         | TCTCGCAG                    | 139           | C06  | ATTGAGCT         | TCACTGTG         | CACAGTGA                    |
| 116           | D03  | ATCATGCG         | CAACCAAC         | GTTGGTTG                    | 140           | D06  | GATAGAGA         | ATTGAGCT         | AGCTCAAT                    |
| 117           | E03  | TCTGAGTC         | ATCATGCG         | CGCATGAT                    | 141           | E06  | TCTAGAGC         | GATAGAGA         | TCTCTATC                    |
| 118           | F03  | TCGCCTGT         | TCTGAGTC         | GACTCAGA                    | 142           | F06  | GAATCGCA         | TCTAGAGC         | GCTCTAGA                    |
| 119           | G03  | GCGCAATT         | TCGCCTGT         | ACAGGCGA                    | 143           | G06  | CTTCACGT         | GAATCGCA         | TGCGATTG                    |
| 120           | H03  | AGACGCCT         | GCGCAATT         | AATTGCGC                    | 144           | H06  | CTCCGTT          | CTTCACGT         | ACGTGAAG                    |

**Table 40** SureSelect Max Index Primer Pairs 145–192, provided in blue 96-well plate

| Primer Pair # | Well | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement | Primer Pair # | Well | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement |
|---------------|------|------------------|------------------|-----------------------------|---------------|------|------------------|------------------|-----------------------------|
| 145           | A07  | TGTGACTA         | CTCCGGTT         | AACCGGAG                    | 169           | A10  | CGCTCAGA         | CTAACAAG         | CTTGTTAG                    |
| 146           | B07  | GCTTCCAG         | TGTGACTA         | TAGTCACA                    | 170           | B10  | TAACGACA         | CGCTCAGA         | TCTGAGCG                    |
| 147           | C07  | CATCCTGT         | GCTTCCAG         | CTGGAAGC                    | 171           | C10  | CATACTTG         | TAACGACA         | TGTCGTTA                    |
| 148           | D07  | GTAATACG         | CATCCTGT         | ACAGGATG                    | 172           | D10  | AGATACGA         | CATACTTG         | CAAGTATG                    |
| 149           | E07  | GCCAACAA         | GTAATACG         | CGTATTAC                    | 173           | E10  | AATCCGAC         | AGATACGA         | TCGTATCT                    |
| 150           | F07  | CATGACAC         | GCCAACAA         | TTGTTGGC                    | 174           | F10  | TGAAGTAC         | AATCCGAC         | GTCGGATT                    |
| 151           | G07  | TGCAATGC         | CATGACAC         | GTGTCATG                    | 175           | G10  | CGAATCAT         | TGAAGTAC         | GTACTIONA                   |
| 152           | H07  | CACATTCG         | TGCAATGC         | GCATTGCA                    | 176           | H10  | TGATTGGC         | CGAATCAT         | ATGATTTCG                   |
| 153           | A08  | CAATCCGA         | CACATTCG         | CGAATGTG                    | 177           | A11  | TCGAAGGA         | TGATTGGC         | GCCAATCA                    |
| 154           | B08  | CATCGACG         | CAATCCGA         | TCGGATTG                    | 178           | B11  | CAGTCATT         | TCGAAGGA         | TCCTTCGA                    |
| 155           | C08  | GTGCGCTT         | CATCGACG         | CGTCGATG                    | 179           | C11  | CGCGAACA         | CAGTCATT         | AATGACTG                    |
| 156           | D08  | ATAGCGTT         | GTGCGCTT         | AAGCGCAC                    | 180           | D11  | TACGGTTG         | CGCGAACA         | TGTTTCGCG                   |
| 157           | E08  | GAGTAAGA         | ATAGCGTT         | AACGCTAT                    | 181           | E11  | AGAACCGT         | TACGGTTG         | CAACCGTA                    |
| 158           | F08  | CTGACACA         | GAGTAAGA         | TCTTACTC                    | 182           | F11  | AGGTGCTT         | AGAACCGT         | ACGGTTCT                    |
| 159           | G08  | ATACGTGT         | CTGACACA         | TGTGTCAG                    | 183           | G11  | ATCGCAAC         | AGGTGCTT         | AAGCACCT                    |
| 160           | H08  | GACCGAGT         | ATACGTGT         | ACACGTAT                    | 184           | H11  | GCCTCTCA         | ATCGCAAC         | GTTGCGAT                    |
| 161           | A09  | GCAGTTAG         | GACCGAGT         | ACTCGGTC                    | 185           | A12  | TCGCGTCA         | GCCTCTCA         | TGAGAGGC                    |
| 162           | B09  | CGTTCGTC         | GCAGTTAG         | CTAACTGC                    | 186           | B12  | GAGTGCGT         | TCGCGTCA         | TGACGCGA                    |
| 163           | C09  | CGTTAACG         | CGTTCGTC         | GACGAACG                    | 187           | C12  | CGAACACT         | GCATAAGT         | ACTTATGC                    |
| 164           | D09  | TCGAGCAT         | CGTTAACG         | CGTTAACG                    | 188           | D12  | TAAGAGTG         | AGAAGACG         | CGTCTTCT                    |
| 165           | E09  | GCCGTAAC         | TCGAGCAT         | ATGCTCGA                    | 189           | E12  | TGGATTGA         | TAAGAGTG         | CACTCTTA                    |
| 166           | F09  | GAGCTGTA         | GCCGTAAC         | GTTACGGC                    | 190           | F12  | AGGACATA         | TGGATTGA         | TCAATCCA                    |
| 167           | G09  | AGGAAGAT         | GAGCTGTA         | TACAGCTC                    | 191           | G12  | GACATCCT         | AGGACATA         | TATGTCCT                    |
| 168           | H09  | CTAACAAG         | AGGAAGAT         | ATCTTCCT                    | 192           | H12  | GAAGCCTC         | GACATCCT         | AGGATGTC                    |

**Table 41** SureSelect Max Index Primer Pairs 193–240, provided in green 96-well plate

| Primer Pair # | Well | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement | Primer Pair # | Well | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement |
|---------------|------|------------------|------------------|-----------------------------|---------------|------|------------------|------------------|-----------------------------|
| 193           | A01  | GTCTCTTC         | GAAGCCTC         | GAGGCTTC                    | 217           | A04  | GCGGTATG         | CACGAGCT         | AGCTCGTG                    |
| 194           | B01  | AGTCACTT         | GTCTCTTC         | GAAGAGAC                    | 218           | B04  | TCTATGCG         | GCGGTATG         | CATACCGC                    |
| 195           | C01  | AGCATACA         | AGTCACTT         | AAGTGACT                    | 219           | C04  | AGGTGAGA         | TCTATGCG         | CGCATAGA                    |
| 196           | D01  | TCAGACAA         | AGCATACA         | TGTATGCT                    | 220           | D04  | CACAACTT         | AGGTGAGA         | TCTCACCT                    |
| 197           | E01  | TTGGAGAA         | TCAGACAA         | TTGTCTGA                    | 221           | E04  | TTGTGTAC         | CACAACTT         | AAGTTGTG                    |
| 198           | F01  | TTAACGTG         | TTGGAGAA         | TTCTCAA                     | 222           | F04  | TCACAAGA         | TTGTGTAC         | GTACACAA                    |
| 199           | G01  | CGTCTGTG         | TTAACGTG         | CACGTAA                     | 223           | G04  | GAAGACCT         | TCACAAGA         | TCTTGTGA                    |
| 200           | H01  | AACCTAAC         | CGTCTGTG         | CACAGACG                    | 224           | H04  | AGTTCTGT         | GAAGACCT         | AGGTCTTC                    |
| 201           | A02  | AGAGTGCT         | AACCTAAC         | GTTAGGTT                    | 225           | A05  | GCAGTGTT         | AGTTCTGT         | ACAGAACT                    |
| 202           | B02  | TTATCTCG         | AGAGTGCT         | AGCACTCT                    | 226           | B05  | AGGCATGC         | GCAGTGTT         | AACACTGC                    |
| 203           | C02  | CATCAGTC         | TTATCTCG         | CGAGATAA                    | 227           | C05  | AAGGTACT         | AGGCATGC         | GCATGCCT                    |
| 204           | D02  | AAGCACAA         | CATCAGTC         | GACTIONG                    | 228           | D05  | CACTAAGT         | AAGGTACT         | AGTACCTT                    |
| 205           | E02  | CAGTGAGC         | AAGCACAA         | TTGTGCTT                    | 229           | E05  | GAGTCCTA         | CACTAAGT         | ACTTAGTG                    |
| 206           | F02  | GTCGAAGT         | CAGTGAGC         | GCTCACTG                    | 230           | F05  | AGTCCTTC         | GAGTCCTA         | TAGGACTC                    |
| 207           | G02  | TCTCATGC         | GTCGAAGT         | ACTTCGAC                    | 231           | G05  | TTAGGAAC         | AGTCCTTC         | GAAGGACT                    |
| 208           | H02  | CAGAAGAA         | TCTCATGC         | GCATGAGA                    | 232           | H05  | AAGTCCAT         | TTAGGAAC         | GTTCCCTAA                   |
| 209           | A03  | CGGATAGT         | CAGAAGAA         | TTCTTCTG                    | 233           | A06  | GAATACGC         | AAGTCCAT         | ATGGACTT                    |
| 210           | B03  | CACGTGAG         | CGGATAGT         | ACTATCCG                    | 234           | B06  | TCCAATCA         | GAATACGC         | GCGTATTC                    |
| 211           | C03  | TACGATAC         | CACGTGAG         | CTCACGTG                    | 235           | C06  | CGACGGTA         | TCCAATCA         | TGATTGGA                    |
| 212           | D03  | CGCATGCT         | TACGATAC         | GTATCGTA                    | 236           | D06  | CATTGCAT         | CGACGGTA         | TACCGTCG                    |
| 213           | E03  | GCTTGCTA         | CGCATGCT         | AGCATGCG                    | 237           | E06  | ATCTGCGT         | CATTGCAT         | ATGCAATG                    |
| 214           | F03  | GAACGCAA         | GCTTGCTA         | TAGCAAGC                    | 238           | F06  | GTACCTTG         | ATCTGCGT         | ACGCAGAT                    |
| 215           | G03  | ATCTACCA         | GAACGCAA         | TTGCGTTC                    | 239           | G06  | GAGCATAC         | GTACCTTG         | CAAGGTAC                    |
| 216           | H03  | CACGAGCT         | ATCTACCA         | TGGTAGAT                    | 240           | H06  | TGCTTACG         | GAGCATAC         | GTATGCTC                    |

**Table 42** SureSelect Max Index Primer Pairs 241–288, provided in green 96-well plate

| Primer Pair # | Well | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement | Primer Pair # | Well | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement |
|---------------|------|------------------|------------------|-----------------------------|---------------|------|------------------|------------------|-----------------------------|
| 241           | A07  | AAGAGACA         | TGCTTACG         | CGTAAGCA                    | 265           | A10  | CAATGCTG         | CATGAATG         | CATTCATG                    |
| 242           | B07  | TAGCTATG         | AAGAGACA         | TGTCTCTT                    | 266           | B10  | CTTGATCA         | CAATGCTG         | CAGCATTG                    |
| 243           | C07  | TCTGCTAC         | TAGCTATG         | CATAGCTA                    | 267           | C10  | GCGAATTA         | CTTGATCA         | TGATCAAG                    |
| 244           | D07  | GTCACAGA         | TCTGCTAC         | GTAGCAGA                    | 268           | D10  | GTTGAGC          | GCGAATTA         | TAATTCGC                    |
| 245           | E07  | CGATTGAA         | GTCACAGA         | TCTGTGAC                    | 269           | E10  | GCCAGTAG         | GTTGAGC          | GCTCGAAC                    |
| 246           | F07  | GAGAGATT         | CGATTGAA         | TTCAATCG                    | 270           | F10  | AAGGTCGA         | GCCAGTAG         | CTACTGGC                    |
| 247           | G07  | TCATACCG         | GAGAGATT         | AATCTCTC                    | 271           | G10  | AGTGAAGT         | CACTTATG         | CATAAGTG                    |
| 248           | H07  | TCCGAACT         | TCATACCG         | CGGTATGA                    | 272           | H10  | GTTGCAAG         | ATAACGGC         | GCCGTTAT                    |
| 249           | A08  | AGAGAGAA         | TCCGAACT         | AGTTCGGA                    | 273           | A11  | AGCCGGAA         | GTTGCAAG         | CTTGCAAC                    |
| 250           | B08  | GATCGTTA         | AGAGAGAA         | TTCTCTCT                    | 274           | B11  | AACAGCCG         | AGCCGGAA         | TTCCGGCT                    |
| 251           | C08  | GCGCTAGA         | GATCGTTA         | TAACGATC                    | 275           | C11  | CTAGTGTA         | AACAGCCG         | CGGCTGTT                    |
| 252           | D08  | ATGACTCG         | GCGCTAGA         | TCTAGCGC                    | 276           | D11  | GAGGCTCT         | CTAGTGTA         | TACTACTAG                   |
| 253           | E08  | CAATAGAC         | ATGACTCG         | CGAGTCAT                    | 277           | E11  | CTCCGCAA         | GAGGCTCT         | AGAGCCTC                    |
| 254           | F08  | CGATATGC         | CAATAGAC         | GTCTATTG                    | 278           | F11  | CGCTATTG         | CTCCGCAA         | TTGCGGAG                    |
| 255           | G08  | GTCAGAAT         | CGATATGC         | GCATATCG                    | 279           | G11  | GTGTTGAG         | CGCTATTG         | CAATAGCG                    |
| 256           | H08  | CATAAGGT         | GCACTACT         | AGTAGTGC                    | 280           | H11  | TCACCGAC         | GTGTTGAG         | CTCAACAC                    |
| 257           | A09  | TGTTGGTT         | GATTCGGC         | GCCGAATC                    | 281           | A12  | CGGTAATC         | TCACCGAC         | GTCGGTGA                    |
| 258           | B09  | ATACTCGC         | TGTTGGTT         | AACCAACA                    | 282           | B12  | GTGACTGC         | CGGTAATC         | GATTACCG                    |
| 259           | C09  | AATGCTAG         | ATACTCGC         | GCGAGTAT                    | 283           | C12  | CGACTTGT         | GTGACTGC         | GCAGTCAC                    |
| 260           | D09  | GCCTAGGA         | AATGCTAG         | CTAGCATT                    | 284           | D12  | GATAGGAC         | CGACTTGT         | ACAAGTCG                    |
| 261           | E09  | GCAACCGA         | GCCTAGGA         | TCCTAGGC                    | 285           | E12  | AAGTACTC         | GATAGGAC         | GTCCTATC                    |
| 262           | F09  | ATACTGCA         | GCAACCGA         | TCGTTGTC                    | 286           | F12  | GCTCTCTC         | AAGTACTC         | GAGTACTT                    |
| 263           | G09  | TCTCCTTG         | ATACTGCA         | TGCAGTAT                    | 287           | G12  | CTACCAGT         | GCTCTCTC         | GAGAGAGC                    |
| 264           | H09  | CATGAATG         | TCTCCTTG         | CAAGGAGA                    | 288           | H12  | GATGAGAT         | CTACCAGT         | ACTGGTAG                    |

**Table 43** SureSelect Max Index Primer Pairs 289–336, provided in red 96-well plate

| Primer Pair # | Well | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement | Primer Pair # | Well | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement |
|---------------|------|------------------|------------------|-----------------------------|---------------|------|------------------|------------------|-----------------------------|
| 289           | A01  | AGATAGTG         | GATGAGAT         | ATCTCATC                    | 313           | A04  | AGCTACAT         | GATCCATG         | CATGGATC                    |
| 290           | B01  | AGAGGTTA         | AGATAGTG         | CACTATCT                    | 314           | B04  | CGCTGTAA         | AGCTACAT         | ATGTAGCT                    |
| 291           | C01  | CTGACCGT         | AGAGGTTA         | TAACCTCT                    | 315           | C04  | CACTACCG         | CGCTGTAA         | TTACAGCG                    |
| 292           | D01  | GCATGGAG         | CTGACCGT         | ACGGTCAG                    | 316           | D04  | GCTCACGA         | CACTACCG         | CGGTAGTG                    |
| 293           | E01  | CTGCCTTA         | GCATGGAG         | CTCCATGC                    | 317           | E04  | TGGCTTAG         | GCTCACGA         | TCGTGAGC                    |
| 294           | F01  | GCGTCACT         | CTGCCTTA         | TAAGGCAG                    | 318           | F04  | TCCAGACG         | TGGCTTAG         | CTAAGCCA                    |
| 295           | G01  | GCGATTAC         | GCGTCACT         | AGTGACGC                    | 319           | G04  | AGTGGCAT         | TCCAGACG         | CGTCTGGA                    |
| 296           | H01  | TCACCACG         | GCGATTAC         | GTAATCGC                    | 320           | H04  | TGTACCGA         | AGTGGCAT         | ATGCCACT                    |
| 297           | A02  | AGACCTGA         | TCACCACG         | CGTGGTGA                    | 321           | A05  | AAGACTAC         | TGTACCGA         | TCGGTACA                    |
| 298           | B02  | GCCGATAT         | AGACCTGA         | TCAGGTCT                    | 322           | B05  | TGCCGTTA         | AAGACTAC         | GTAGTCTT                    |
| 299           | C02  | CTTATTGC         | GCCGATAT         | ATATCGGC                    | 323           | C05  | TTGGATCT         | TGCCGTTA         | TAACGGCA                    |
| 300           | D02  | CGATACCT         | CTTATTGC         | GCAATAAG                    | 324           | D05  | TCCTCAA          | TTGGATCT         | AGATCCAA                    |
| 301           | E02  | CTCGACAT         | CGATACCT         | AGGTATCG                    | 325           | E05  | CGAGTCGA         | TCCTCAA          | TTGGAGGA                    |
| 302           | F02  | GAGATCGC         | CTCGACAT         | ATGTGAG                     | 326           | F05  | AGGCTCAT         | CGAGTCGA         | TCGACTCG                    |
| 303           | G02  | CGGTCTCT         | GAGATCGC         | GCGATCTC                    | 327           | G05  | GACGTGCA         | AGGCTCAT         | ATGAGCCT                    |
| 304           | H02  | TAACAC           | CGGTCTCT         | AGAGACCG                    | 328           | H05  | GAACATGT         | GACGTGCA         | TGCACGTC                    |
| 305           | A03  | CACAATGA         | TAACAC           | GTGAGTTA                    | 329           | A06  | AATTGGCA         | GAACATGT         | ACATGTTC                    |
| 306           | B03  | GACTGACG         | CACAATGA         | TCATTGTG                    | 330           | B06  | TGGAGACT         | AATTGGCA         | TGCCAATT                    |
| 307           | C03  | CTTAAGAC         | GACTGACG         | CGTCAGTC                    | 331           | C06  | AACTCACA         | TGGAGACT         | AGTCTCCA                    |
| 308           | D03  | GAGTGTAG         | CTTAAGAC         | GTCTTAAG                    | 332           | D06  | GTAGACTG         | AACTCACA         | TGTGAGTT                    |
| 309           | E03  | TGCACATC         | GAGTGTAG         | CTACACTC                    | 333           | E06  | CGTAGTTA         | GTAGACTG         | CAGTCTAC                    |
| 310           | F03  | CGATGTCG         | TGCACATC         | GATGTGCA                    | 334           | F06  | CGTCAGAT         | CGTAGTTA         | TAACACG                     |
| 311           | G03  | AACACCGA         | CGATGTCG         | CGACATCG                    | 335           | G06  | AACGGTCA         | CGTCAGAT         | ATCTGACG                    |
| 312           | H03  | GATCCATG         | AACACCGA         | TCGGTGTT                    | 336           | H06  | GCCTTCAT         | AACGGTCA         | TGACCGTT                    |

**Table 44** SureSelect Max Index Primer Pairs 337–384, provided in red 96-well plate

| Primer Pair # | Well | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement | Primer Pair # | Well | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement |
|---------------|------|------------------|------------------|-----------------------------|---------------|------|------------------|------------------|-----------------------------|
| 337           | A07  | TGAGACGC         | GCCTTCAT         | ATGAAGGC                    | 361           | A10  | CTGAGCTA         | GCACAGTA         | TACTGTGC                    |
| 338           | B07  | CATCGGAA         | TGAGACGC         | GCGTCTCA                    | 362           | B10  | CTTGCGAT         | CTGAGCTA         | TAGCTCAG                    |
| 339           | C07  | TAGGACAT         | CATCGGAA         | TTCCGATG                    | 363           | C10  | GAAGTAGT         | CTTGCGAT         | ATCGCAAG                    |
| 340           | D07  | AACACAAG         | TAGGACAT         | ATGTCCTA                    | 364           | D10  | GTTATCGA         | GAAGTAGT         | ACTACTTC                    |
| 341           | E07  | TTCGACTC         | AACACAAG         | CTTGTGTT                    | 365           | E10  | TGTCGTCG         | GTTATCGA         | TCGATAAC                    |
| 342           | F07  | GTCGGTAA         | TTCGACTC         | GAGTCGAA                    | 366           | F10  | CGTAACTG         | TGTCGTCG         | CGACGACA                    |
| 343           | G07  | GTTTCATTC        | GTCGGTAA         | TTACCGAC                    | 367           | G10  | GCATGCCT         | CGTAACTG         | CAGTTACG                    |
| 344           | H07  | AAGCAGTT         | GTTTCATTC        | GAATGAAC                    | 368           | H10  | TCGTACAC         | GCATGCCT         | AGGCATGC                    |
| 345           | A08  | ATAAGCTG         | AAGCAGTT         | AACTGCTT                    | 369           | A11  | CACAGGTG         | TCGTACAC         | GTGTACGA                    |
| 346           | B08  | GCTTAGCG         | ATAAGCTG         | CAGCTTAT                    | 370           | B11  | AGCAGTGA         | CACAGGTG         | CACCTGTG                    |
| 347           | C08  | TTCCAACA         | GCTTAGCG         | CGCTAAGC                    | 371           | C11  | ATTCCAGA         | AGCAGTGA         | TCACTGCT                    |
| 348           | D08  | TACCGCAT         | TTCCAACA         | TGTTGGAA                    | 372           | D11  | TCCTTGAG         | ATTCCAGA         | TCTGGAAT                    |
| 349           | E08  | AGGCAATG         | TACCGCAT         | ATGCGGTA                    | 373           | E11  | ATACCTAC         | TCCTTGAG         | CTCAAGGA                    |
| 350           | F08  | GCCTCGTT         | AGGCAATG         | CATTGCCT                    | 374           | F11  | AGACCATT         | ATACCTAC         | GTAGGTAT                    |
| 351           | G08  | CACGGATC         | GCCTCGTT         | AACGAGGC                    | 375           | G11  | CGTAAGCA         | AGACCATT         | AATGGTCT                    |
| 352           | H08  | GAGACACG         | CACGGATC         | GATCCGTG                    | 376           | H11  | TCTGTCAG         | CGTAAGCA         | TGCTTACG                    |
| 353           | A09  | AGAGTAAG         | GAGACACG         | CGTGTCTC                    | 377           | A12  | CACAGACT         | TCTGTCAG         | CTGACAGA                    |
| 354           | B09  | AGTACGTT         | AGAGTAAG         | CTTACTCT                    | 378           | B12  | GTCGCCTA         | CACAGACT         | AGTCTGTG                    |
| 355           | C09  | AACGCTGC         | AGTACGTT         | AACGTAAT                    | 379           | C12  | TGCGCTCT         | GTCGCCTA         | TAGGCGAC                    |
| 356           | D09  | GTAGAGCA         | AACGCTGC         | GCAGCGTT                    | 380           | D12  | GCTATAAG         | TGCGCTCT         | AGAGCGCA                    |
| 357           | E09  | TCCTGAGA         | GTAGAGCA         | TGCTCTAC                    | 381           | E12  | CAACAAC          | GCTATAAG         | CTTATAGC                    |
| 358           | F09  | CTGAATAG         | TCCTGAGA         | TCTCAGGA                    | 382           | F12  | AGAGAATC         | CTCTCACT         | AGTGAGAG                    |
| 359           | G09  | CAAGACTA         | CTGAATAG         | CTATTCAG                    | 383           | G12  | TAATGGTC         | AGACGAGC         | GCTCGTCT                    |
| 360           | H09  | GCACAGTA         | CAAGACTA         | TAGTCTTG                    | 384           | H12  | GTTGTATC         | TAATGGTC         | GACCATTA                    |

# 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 yield of libraries is low

- ✓ Use only nuclease-free water or 1X Low TE Buffer (10 mM Tris-HCl pH 8.0, 0.1 mM EDTA) where specified in the protocol for sample preparation and dilution steps. Do not substitute with standard TE or other solvents containing >0.1 mM EDTA. Higher concentrations of EDTA inhibit the fragmentation reaction and reduce library yield.
- ✓ 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.
- ✓ PCR cycle number may require optimization. Repeat library preparation for the sample, increasing the 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 over-amplified. Repeat library preparation for the sample, decreasing the PCR cycle number by 1 to 3 cycles.
- ✓ DNA isolated from 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.
- ✓ Yield from the magnetic bead purification step may be suboptimal. Consider the factors below:
  - Adhere to all of the bead and reagent handling steps in the protocol. In particular, make sure to equilibrate beads at room temperature for at least 30 minutes before use, and use freshly-prepared 70% ethanol (prepared on day of use) for each purification procedure.
  - Ensure that the magnetic beads are not over-dried just prior to sample elution (see [step 12 on page 20](#) and [page 23](#)). Monitor the bead pellets during the drying incubation frequently and conclude the drying step immediately after the residual ethanol has evaporated.
  - Increasing the elution incubation time (up to 10 minutes) may improve recovery, especially for longer DNA fragments.

## If library fragment size is different than expected or required in electropherograms

- ✓ If library fragments are longer than expected the gDNA samples may be under-fragmented. Use only nuclease-free water or 1X Low TE Buffer (10 mM Tris-HCl pH 8.0, 0.1 mM EDTA) where specified in the protocol for sample preparation and dilution steps. Do not substitute with standard TE or other solvents containing >0.1 mM EDTA. Higher concentrations of EDTA can inhibit DNA fragmentation.
- ✓ If library fragments are shorter than expected the gDNA samples may be over-fragmented. Make sure to keep DNA samples on ice while setting up the enzymatic fragmentation/end-repair/dA-tailing reactions in [step 3 on page 18](#).
- ✓ FFPE DNA pre-capture libraries may have a smaller fragment size distribution due to the presence of DNA fragments in the sample input that are smaller than the targeted post-fragmentation size. Adhere to the DNA quality guidelines provided on [page 16](#).

- ✓ DNA fragment size selection during magnetic bead purification depends upon using the correct ratio of sample to purification beads. Before removing an aliquot of beads for the purification step, mix the beads until the suspension appears homogeneous and consistent in color. Verify that you are using the recommended bead volume at each stage of the protocol.
- ✓ For the whole genome sequencing workflow, the average library fragment size can be increased by adjusting the ratio of bead suspension to DNA sample volumes in the final library purification procedure on [page 30](#) as detailed below. Note that this protocol modification will also decrease the overall library yield.
  - For 2x150 NGS the standard protocol uses beads at 1X the library DNA sample volume. To increase the average fragment size in the library, decrease the bead volume ratio to the range of 0.6X to 0.7X (with 30  $\mu$ L or 35  $\mu$ L of purification bead suspension added to each 50  $\mu$ L sample).
  - For 2x250 NGS the standard protocol uses beads at 0.6X the library DNA sample volume. To increase the average fragment size in the library, decrease the bead volume ratio to the range of 0.5X to 0.55X (with 50  $\mu$ L or 55  $\mu$ L of purification bead suspension added to each 100  $\mu$ L sample).

### **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 25](#). 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 19](#). In particular, ensure that the Adaptor Oligo Mix is mixed with the sample prior to adding the Ligation Master Mix to the mixture. Do not add the Ligation Master Mix and the Adaptor Oligo Mix to the sample in a single step.

## Quick Reference Protocol: Library Prep for Target Enrichment

An abbreviated protocol summary is provided below for experienced users. Use the complete protocol on [page 14](#) to [page 25](#) until you are familiar with all of the protocol details.

| Step   | Summary of Conditions  |
|--|--|
| Prepare and qualify DNA samples                  | Prepare 10–200 ng gDNA in 40 µL nuclease-free H <sub>2</sub> O or 1X Low TE Buffer> place in sample wells> keep on ice.<br>For FFPE DNA, qualify integrity and adjust input amount as directed on <a href="#">page 16</a> .  |
| Prepare Frag/A-Tail master mix                   | <b>Per 8 reactions:</b> 36 µL Frag/A-Tail Buffer + 54 µL Frag/A-Tail Enzyme Mix<br><b>Per 24 reactions:</b> 104 µL Frag/A-Tail Buffer + 156 µL Frag/A-Tail Enzyme Mix<br>Prepare on ice> vortex> spin> keep on ice.  |
| Fragment, end-repair and dA-Tail the DNA samples | 40 µL DNA sample + 10 µL Frag/A-Tail master mix<br>Mix> spin> incubate in thermal cycler using program in <a href="#">Table 45</a> .   |
| Ligate adaptor                                   | 50 µL DNA fragments + 5 µL SureSelect Max Adaptor Oligo Mix (MBC or MBC-free)> mix> spin.<br>Add 20 µL Ligation Master Mix> pipette to mix> spin.<br>Incubate in thermal cycler: Hold @ 4°C until samples loaded, 15 min @ 20°C, Hold @ 4°C.   |
| Purify DNA                                       | 75 µL DNA sample + 60 µL purification bead suspension> mix> incubate 5 min> collect beads> discard supernatant. Wash beads 2X with 70% ethanol.<br>Elute DNA in 21 µL nuclease-free H <sub>2</sub> O> mix> incubate 2-5 minutes> collect beads> transfer 20 µL supernatant to fresh well.    |
| Index and amplify library                        | 20 µL DNA sample + 25 µL Amplification Master Mix + 5 µL SureSelect Max UDI Primers for ILM<br>Vortex> spin> amplify in thermal cycler using program in <a href="#">Table 46</a> .   |
| Purify amplified library DNA                     | 50 µL amplified DNA + 50 µL purification bead suspension> mix> incubate 5 min> collect beads> discard supernatant. Wash beads 2X with 70% ethanol.<br>Elute DNA in 15 µL nuclease-free H <sub>2</sub> O> mix> incubate 2-5 minutes> collect beads> transfer 14 µL supernatant to fresh well. |
| Quantify and qualify DNA                         | Optional: Analyze quantity and quality using TapeStation or Fragment Analyzer System   |

**Table 45** Frag/A-Tail reaction thermal cycler program for target enrichment workflow (50 µL vol, heated lid at 105°C)

| Segment | Temperature | Time   |
|---------|-------------|--|
| 1       | 4°C         | Hold—proceed to Step 2 once samples added to block   |
| 2       | 37°C        | Intact DNA for 2 ×100 NGS: 20 minutes<br>Intact DNA for 2 ×150 NGS: 10 minutes<br>FFPE DNA for any read length NGS: 15 minutes |
| 3       | 65°C        | 30 minutes   |
| 4       | 4°C         | Hold   |

**Table 46** Library indexing/amplification thermal cycler program for target enrichment workflow (50 µL vol; heated lid at 105°C)

| Segment | Number of Cycles  | Temperature   | Time   |
|---------|---|---|--|
| 1       | 1   | 98°C  | 45 seconds   |
| 2       | Intact DNA input<br>200 ng: 7 cycles<br>100 ng: 8 cycles<br>50 ng: 9 cycles<br>10 ng: 10 cycles | OR<br>FFPE DNA input<br>100–200 ng: 11 cycles<br>50 ng: 12 cycles<br>10 ng: 13 cycles | 98°C<br>15 seconds<br>60°C<br>30 seconds<br>72°C<br>30 seconds |
| 3       | 1   | 72°C  | 1 minute   |
| 4       | 1   | 4°C   | Hold   |

# Quick Reference Protocol: Library Prep for Whole Genome NGS

An abbreviated protocol summary is provided below for experienced users. Use the complete protocol on [page 28](#) to [page 32](#) until you are familiar with all of the protocol details.

| Step   | Summary of Conditions  |
|--|--|
| Prepare and qualify DNA samples                  | Prepare 10–200 ng gDNA in 40 µL 1X Low TE Buffer (recommended) or nuclease-free H <sub>2</sub> O> place in sample wells> keep on ice.<br>For FFPE DNA, qualify integrity and adjust input amount as directed on <a href="#">page 16</a> .  |
| Prepare Frag/A-Tail master mix                   | <b>Per 8 reactions:</b> 36 µL Frag/A-Tail Buffer + 54 µL Frag/A-Tail Enzyme Mix<br><b>Per 24 reactions:</b> 104 µL Frag/A-Tail Buffer + 156 µL Frag/A-Tail Enzyme Mix<br>Prepare on ice> vortex> spin> keep on ice.  |
| Fragment, end-repair and dA-Tail the DNA samples | 40 µL DNA sample + 10 µL Frag/A-Tail master mix<br>Mix> spin> incubate in thermal cycler using program in <a href="#">Table 47</a> .   |
| Ligate adaptor                                   | 50 µL DNA fragments + 5 µL SureSelect Max MBC-Free Adaptor Oligo Mix> mix> spin.<br>Add 20 µL Ligation Master Mix> pipette to mix> spin.<br>Incubate in thermal cycler: Hold @ 4°C until samples loaded, 15 min @ 20°C, Hold @ 4°C.  |
| Purify DNA                                       | 75 µL DNA sample + 60 µL purification bead suspension> mix> incubate 5 min> collect beads> discard supernatant. Wash beads 2X with 70% ethanol.<br>Elute DNA in 21 µL nuclease-free H <sub>2</sub> O> mix> incubate 2-5 minutes> collect beads> transfer 20 µL supernatant to fresh well.  |
| Index and amplify library                        | 20 µL DNA sample + 25 µL Amplification Master Mix + 5 µL SureSelect Max UDI Primers for ILM<br>Vortex> spin> amplify in thermal cycler using program in <a href="#">Table 48</a> .   |
| For 2x250 only, dilute 2X                        | For libraries to be sequenced using 2x250 reads, add 50 µL of nuclease-free H <sub>2</sub> O per sample well.  |
| Purify amplified library DNA                     | <b>For 2x150 samples:</b> 50 µL amplified DNA + 50 µL purification bead suspension<br><b>For 2x250 samples:</b> 100 µL amplified DNA dilution + 60 µL purification bead suspension<br>Mix> incubate 10 min> collect beads> discard supernatant. Wash beads 2X with 70% ethanol.<br>Elute DNA in 26 µL Low TE Buffer> mix> incubate 2-5 minutes> collect beads> transfer 25 µL supernatant to fresh well. |
| Quantify and qualify DNA                         | QC by TapeStation or Fragment Analyzer. Quantify library DNA for NGS using qPCR for 3- or 4-cycle amplified libraries or using TapeStation or Fragment Analyzer for libraries amplified with ≥5 cycles.  |

**Table 47** Frag/A-Tail reaction thermal cycler program for whole genome workflow (50 µL vol, heated lid at 105°C)

| Segment | Temperature | Time  |
|---------|-------------|---|
| 1       | 4°C         | Hold--proceed to Segment 2 once samples added to block  |
| 2       | 37°C        | Intact DNA for 2x150 NGS: 5 minutes<br>Intact DNA for 2x250 NGS: 2 minutes<br>FFPE DNA for any read length NGS: 5 minutes |
| 3       | 65°C        | 30 minutes  |
| 4       | 4°C         | Hold  |

**Table 48** Library indexing/amplification thermal cycler program for whole genome workflow (50  $\mu$ L vol; heated lid at 105°C)

| Segment | Number of Cycles   | Temperature | Time       |
|---------|--|-------------|------------|
| 1       | 1  | 98°C        | 45 seconds |
| 2       | 50–200 ng Intact DNA input libraries: 3–5 cycles<br>10 ng Intact DNA input libraries: 6 cycles | 98°C        | 15 seconds |
|         |  | 60°C        | 30 seconds |
|         | 50–200 ng FFPE DNA input libraries: 7 cycles<br>10 ng FFPE DNA input libraries: 8 cycles       | 72°C        | 30 seconds |
| 3       | 1  | 72°C        | 1 minute   |
| 4       | 1  | 4°C         | Hold       |

## In This Book

This guide provides instructions for SureSelect Max DNA Library Preparation with Enzymatic Fragmentation for Illumina sequencing.

