



SureSelect Max Target Enrichment using Overnight Hybridization

For NGS using the Illumina Platform

Protocol

Version A0 September 2024

SureSelect platform manufactured with Agilent SurePrint technology.

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

Notices

© Agilent Technologies, Inc. 2024

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

G9690-90000

Edition

Version A0 September 2024

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

Acknowledgment

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

Technical Support

For US and Canada

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

Or send an e-mail to:

ngs.support@agilent.com

For all other regions

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

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

Warranty

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

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 target enrichment of Illumina paired-end multiplexed DNA or cDNA sequencing libraries using SureSelect Max Overnight Hybridization Modules and a SureSelect probe. Libraries can be pooled for NGS using either pre-capture or post-capture pooling. The target-enriched libraries are ready for NGS using a suitable Illumina instrument. Before starting this protocol, SureSelect Max DNA or RNA libraries must be prepared as described in separate guides for earlier workflow segments.

1 Before You Begin

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

2 Target Enrichment with Overnight Hybridization Protocol

This section describes the steps to hybridize and capture targeted fragments in a SureSelect Max library using a SureSelect Probe, using overnight (16–24 hour) hybridization conditions.

3 Appendix: NGS and Analysis Guidelines

This section provides guidelines for downstream NGS and analysis, including information on the SureSelect Max UDIs added to fragments during library preparation.

4 Reference

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

Content

| | | |
|----------|--|-----------|
| 1 | Before You Begin | 5 |
| | Overview of the Workflow | 6 |
| | SureSelect Max Modules and Probes Used in the Workflow | 7 |
| | Additional Materials Used in the Workflow | 8 |
| | Procedural and Safety Notes | 9 |
| 2 | Target Enrichment with Overnight Hybridization Protocol | 10 |
| | Step 1. Place libraries or library pools in the hybridization wells | 11 |
| | Option 1: Plate individual libraries (post-capture pooling workflow) | 11 |
| | Option 2: Plate library pools (pre-capture pooling workflow) | 12 |
| | Step 2. Hybridize libraries to the SureSelect probe | 13 |
| | Step 3. Prepare streptavidin beads and buffers for capture | 16 |
| | Step 4. Capture the hybridized libraries | 16 |
| | Step 5. Amplify the captured libraries | 18 |
| | Step 6. Purify the final libraries using magnetic purification beads | 20 |
| | Step 7. QC and quantify final libraries | 21 |
| | Post-capture pooling guidelines | 23 |
| 3 | Appendix: NGS and Analysis Guidelines | 24 |
| | SureSelect Max Library Composition | 25 |
| | Sequencing Setup and Run Guidelines | 26 |
| | Analysis Pipeline Guidelines | 27 |
| | AGeNT software guidelines for SureSelect Max DNA or RNA workflows | 28 |
| | RNA strandedness guidelines | 28 |
| | SureSelect Max UDI Information | 28 |
| | SureSelect Max index sequences | 30 |
| 4 | Reference | 38 |
| | Reagent Kit Contents | 39 |
| | Component Kit Details | 39 |
| | Troubleshooting Guide | 40 |
| | Quick Reference Protocol: Max Overnight Hyb Target Enrichment | 42 |

1

Before You Begin

Overview of the Workflow [6](#)

SureSelect Max Modules and Probes Used in the Workflow [7](#)

Additional Materials Used in the Workflow [8](#)

Procedural and Safety Notes [9](#)

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 target enrichment workflow segment, using overnight hybridization conditions for target enrichment of prepared SureSelect Max DNA or RNA libraries with a compatible SureSelect XT probe, summarized in [Figure 1](#). The overnight hybridization protocol includes options for post-capture or pre-capture library pooling.

For detailed protocols see [“Target Enrichment with Overnight Hybridization Protocol”](#) on page 10. Protocols for the upstream DNA or RNA library preparation modules are provided in separate publications.

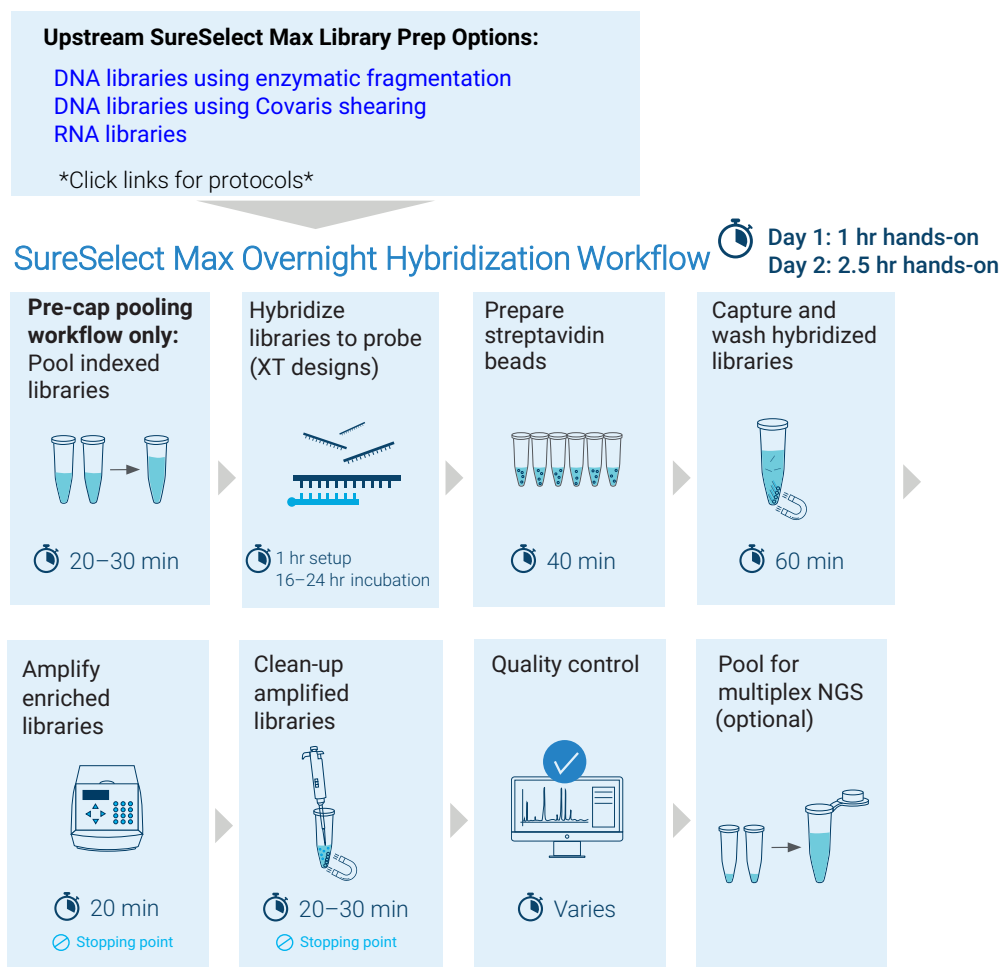


Figure 1 Summary of the SureSelect Max Overnight Hyb target enrichment workflow segment, including optional stopping points and estimated time requirements. Estimates are guidelines for 16 reaction runs. Timing for runs using different protocol parameters may vary.

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

SureSelect Max Modules and Probes Used in the Workflow

This publication provides optimized protocols for the target enrichment workflow segment, using overnight hybridization conditions for target enrichment of prepared SureSelect Max DNA or RNA libraries with a SureSelect XT probe. Agilent's SureSelect reagents required to complete the protocols are summarized in [Table 1](#) and [Table 2](#).

Table 1 SureSelect Max kits used in the Target Enrichment with Overnight Hybridization workflow

| Kit Description | 16 Hybridization Kits* | 96 Hybridization Kits |
|---|---|-----------------------|
| SureSelect Max Overnight Hyb Kit [†] | G9690A | G9690B |
| SureSelect Max Blockers and Primers Kit for ILM | G9699A | G9699B |
| SureSelect Max Purification Beads [‡] | G9962A (5 mL) | G9962B (30 mL) |
| SureSelect XT Probe | See Table 2 for ordering information. Select Overnight Hyb/XT design probes for use in the SureSelect Max Overnight Hybridization workflow. Select the appropriate formulation and format for your indexed library pooling workflow choice (post-capture or pre-capture pooling). | |

* 16 Hyb Kits can be used to process 16 prepared libraries in a post-capture pooling workflow, or up to 96 prepared libraries in a pre-capture pooling workflow.

† Includes SureSelect Streptavidin Beads. Separate purchase not required.

‡ May be substituted with AMPure XP beads (see [Table 3](#) on page 8).

Table 2 Recommended Probes

| Probe Description | Design ID | Ordering Information |
|--|-----------|--|
| Pre-designed Probes | | |
| SureSelect XT Human All Exon V8 | S33266436 | Please visit the Agilent.com probe webpages or the SureDesign website to obtain ordering information for pre-designed probes formulated for either post-capture pooling or pre-capture pooling. Please contact Sales or your local representative if you need assistance. |
| SureSelect XT Human All Exon V8+UTR | S33613367 | |
| SureSelect XT Human All Exon V8+NCV | S33700246 | |
| SureSelect XT Human All Exon V7 | S31285117 | |
| SureSelect XT Clinical Research Exome V4 | S34226363 | |
| SureSelect XT Clinical Research Exome V2 | S30409818 | |
| Custom Probes | | |
| SureSelect Custom Tier1 1–499 kb | | Please visit the SureDesign website to design Custom SureSelect probes and obtain ordering information for custom probes formulated for either post-capture pooling or pre-capture pooling. Contact the SureSelect support team (see page 2) or your local representative if you need assistance. |
| SureSelect Custom Tier2 0.5 –2.9 Mb | | |
| SureSelect Custom Tier3 3 –5.9 Mb | | |
| SureSelect Custom Tier4 6 –11.9 Mb | | |
| SureSelect Custom Tier5 12–24 Mb | | |
| Agilent Community Designs: Please visit the Community Designs (NGS) webpages at agilent.com for information on custom panels developed in collaboration with experts in various fields. | | Design details and ordering information are available at the SureDesign website on the <i>Published Designs</i> tab. Contact the SureSelect support team (see page 2) or your local representative if you need assistance. |

Additional Materials Used in the Workflow

See [Table 3](#) and [Table 4](#) for additional reagents and equipment used in the workflow.

CAUTION

Sample volumes exceed 0.2 mL in certain steps of this protocol. Make sure that the plasticware used with the selected thermal cycler holds ≥ 0.25 mL per well (see [Table 3](#)).

Table 3 Ordering Information for Additional Reagents and Equipment

| Description | Vendor and Part Number | Usage Notes |
|--|---|--|
| Thermal Cyclers (2) with 96-well, 0.2 mL blocks | Various suppliers | Protocols require two thermal cyclers to complete a reagent pre-warming step during certain sample incubation steps. See <i>Note</i> on page 13 for more information. |
| Plasticware compatible with the selected thermal cycler: 8-well strip tubes Tube cap strips (flat or domed)* | Consult the thermal cycler manufacturer's recommendations | Runs using 1–3 strip tubes per run are recommended to facilitate efficient sample handling and to minimize the number of required thermal cyclers. Runs may also be completed using 96 well plates. |
| Vacuum concentrator | Savant SpeedVac, model DNA120, or equivalent | Used for sample dehydration prior to hybridization (4 μ L sample volume required). If a vacuum concentrator is not available in your laboratory, see <i>Troubleshooting</i> on page 40 for suggested protocol modifications. |
| 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 device configured to collect the particles in ring formation. |
| 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 | — |
| Nucleic acid analysis system | See Table 4 on page 9 | — |
| 1X Low TE Buffer | Thermo Fisher Scientific p/n 12090-015, or equivalent | 10 mM Tris-HCl, pH 7.5-8.0, 0.1 mM EDTA |
| 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: AMPure XP Kit (5 mL) | Beckman Coulter Genomics p/n A63880 | Optional alternative to SureSelect Max Purification Beads (See Table 1) |

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

Table 4 Recommended Nucleic Acid Analysis Platforms*

| Analysis System | Vendor and Part Number Information |
|---|-------------------------------------|
| Agilent 4200/4150 TapeStation Instrument | Agilent p/n G2991AA/G2992AA |
| Consumables: | |
| 8-well tube strips | p/n 401428 |
| 8-well tube strip caps | p/n 401425 |
| High Sensitivity D1000 ScreenTape | p/n 5067-5584 |
| High Sensitivity D1000 Reagents | p/n 5067-5585 |
| Agilent 5200/5300/5400 Fragment Analyzer Instrument | Agilent p/n M5310AA/M5311AA/M5312AA |
| Consumables: | |
| HS NGS Fragment Kit (1-6000 bp) | p/n DNF-474-0500 |

* Libraries may also be analyzed using the Agilent 2100 Bioanalyzer instrument, p/n G2939BA, and the High Sensitivity DNA Kit (p/n 5067-4626).

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.
- 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 Target Enrichment with Overnight Hybridization Protocol

- Step 1. Place libraries or library pools in the hybridization wells 11
- Step 2. Hybridize libraries to the SureSelect probe 13
- Step 3. Prepare streptavidin beads and buffers for capture 16
- Step 4. Capture the hybridized libraries 16
- Step 5. Amplify the captured libraries 18
- Step 6. Purify the final libraries using magnetic purification beads 20
- Step 7. QC and quantify final libraries 21
- Post-capture pooling guidelines 23

In this workflow segment, SureSelect Max DNA or cDNA libraries are hybridized with a target-specific probe. After hybridization, the targeted molecules are captured on streptavidin beads and then PCR-amplified. The libraries can be pooled for multiplex NGS after hybridization (see [page 11](#) for post-capture pooling workflow setup) or prior to hybridization (see [page 12](#) for pre-capture pooling workflow setup).

CAUTION

During the overnight hybridization incubation, wells must be adequately sealed to minimize evaporation from the sample wells, or your results can be negatively impacted.

Before you do the first experiment, make sure the plasticware and capping method are appropriate for the thermal cycler. Check that no more than 4 μ L is lost to evaporation under the conditions used for hybridization (see [page 15](#)).

The hybridization workflow segment uses the components listed in [Table 5](#). Prepare as directed before use (refer to the *Where Used* column).

Table 5 Reagents for Hybridization and Capture

| Storage Location | Kit Component | Preparative Steps | Where Used |
|--|---|---|-------------------------|
| -80°C | SureSelect Probe | Thaw on ice before starting Hyb setup on page 13 and keep on ice; vortex to mix | page 15 |
| SureSelect Max Target Enrichment Kit Overnight Hyb Module Box 1, stored at RT | SureSelect Hyb 1 (orange cap or bottle) | Keep at RT, vortex to mix | page 13 |
| | SureSelect Hyb 2 (red cap) | Keep at RT, vortex to mix | page 13 |
| SureSelect Max Target Enrichment Kit Overnight Hyb Module Box 2, stored at -20°C | SureSelect Hyb 4 (black cap) | Keep at RT, vortex to mix | page 13 |
| | SureSelect Hyb 3 (yellow cap) | Thaw and keep at RT, vortex to mix | page 13 |
| SureSelect Max Blockers and Primers Module for ILM, stored at -20°C | SureSelect RNase Block (purple cap) | Thaw and keep on ice, vortex to mix | page 14 |
| | Blocker Mix, ILM (blue cap) | Thaw and keep on ice, vortex to mix | page 14 |

Step 1. Place libraries or library pools in the hybridization wells

Option 1: Plate individual libraries (post-capture pooling workflow)

The hybridization reaction requires prepared library samples in 4 µL of nuclease-free water. See [Table 6](#) for the recommended and required amounts of prepared library sample input for hybridization. To maximize complexity of the final enriched NGS library, add the full 12 µL of prepared library sample to the hybridization reaction. After plating, sample volumes are reduced to 4 µL as described below.

Agilent recommends retaining the remaining prepared library solution volume (~2 µL) for any later troubleshooting QC needs.

Table 6 Hybridization input guidelines

| Type of prepared library | Recommended amount | Minimum amount* |
|--------------------------|---|----------------------|
| gDNA library (DNA input) | 12 µL of undiluted prepared gDNA library solution (pre-capture QC optional) reduced to 4 µL | ≥500 ng library DNA |
| cDNA library (RNA input) | 12 µL of undiluted prepared cDNA library solution (pre-capture QC optional) reduced to 4 µL | ≥200 ng library cDNA |

* For workflows that include the optional pre-capture QC step, use at least the minimum amount of library shown here in the hybridization reactions. When library normalization in the run is not required, using the maximum amount of each library (12 µL of undiluted sample) is still recommended to maximize complexity.

- 1 Prepare the hybridization strip(s) by placing the appropriate amount of each prepared library into a separate sample well.
- 2 Using a vacuum concentrator at ≤45°C, dehydrate the library samples to volume <4 µL without completely drying the samples. Measure the final volume in each well and then bring the volume to 4 µL with nuclease-free water.

NOTE

If a vacuum concentrator is not available in your laboratory, see *Troubleshooting* on [page 40](#).

- 3 Cap the wells, then vortex the strip vigorously for 5–10 seconds. Spin briefly to collect the liquid, then keep the samples on ice until use in [step 4](#) on [page 14](#).

Option 2: Plate library pools (pre-capture pooling workflow)

- 1 Pool the indexed library samples as directed in [Table 7](#), using equal amounts of 8 or 16 libraries, based on the probe design. For each pool, combine the appropriate volume of each prepared library sample in one well of a strip tube.

Table 7 Pre-capture pooling recommendations

| Probe description | Number of indexes per pool | Library type | Total amount of library DNA per pool | Amount of each library |
|--|----------------------------|------------------|--------------------------------------|------------------------|
| SureSelect XT PreCap Human All Exon and Exome probes | 8 | DNA | 4 µg | 500 ng |
| | 8 | cDNA (RNA input) | 1.6 µg | 200 ng |
| SureSelect XT PreCap Custom Probes | 16 | DNA | 4 µg | 250 ng |
| | 16 | cDNA (RNA input) | 1.6 µg | 100 ng |

- 2 Using a vacuum concentrator at $\leq 45^{\circ}\text{C}$, dehydrate the prepped library pools to volume $< 4\ \mu\text{L}$ without completely drying the samples. Measure the final volume in each well and then bring the volume of each pool to $4\ \mu\text{L}$ with nuclease-free water.

NOTE

If a vacuum concentrator is not available in your laboratory, see *Troubleshooting* on [page 40](#).

- 3 Cap the wells, then vortex the sample strip vigorously for 5–10 seconds. Spin briefly to collect the liquid, then keep the samples on ice until use in [step 4](#) on [page 14](#).

Step 2. Hybridize libraries to the SureSelect probe

NOTE

Two thermal cyclers are required for the Hybridization and Capture protocols in the following sections, to complete reagent pre-warming steps during two sample incubation steps: 1) capture bead pre-warming in [step 6](#) on [page 16](#) during hybridization and 2) Wash Buffer 2 pre-warming during the capture incubation (see [step 2](#) and [step 3](#) on [page 17](#)). If two thermal cyclers are not available in your laboratory, the reagent pre-warming steps may be completed using an alternative temperature control device. Reaction mixtures that include the library samples should be incubated using a thermal cycler where directed.

- 1 Prepare the Hybridization Buffer by mixing the components in [Table 8](#) at room temperature. Keep at room temperature until use in [step 7](#). If a precipitate forms, warm the Hybridization Buffer at 65°C for 5 minutes.

Table 8 Preparation of Hybridization Buffer

| Reagent | Volume for 1 reaction | Volume for 8 Hyb reactions (includes excess)* | Volume for 24 Hyb reactions (includes excess)† |
|--------------------------------------|-----------------------|---|--|
| SureSelect Hyb 1 (orange cap/bottle) | 6.63 µL | 66.3 µL | 185.6 µL |
| SureSelect Hyb 2 (red cap) | 0.27 µL | 2.7 µL | 7.6 µL |
| SureSelect Hyb 3 (yellow cap) | 2.65 µL | 26.5 µL | 74.2 µL |
| SureSelect Hyb 4 (black cap) | 3.45 µL | 34.5 µL | 96.6 µL |
| Total | 13 µL | 130 µL | 364 µL |

* 16-Hyb Target Enrichment Kits contain enough reagents for 2 runs of 8 Hybs each using the indicated excess volume.

† 96-Hyb Target Enrichment Kits contain enough reagents for 4 runs of 24 Hybs each using the indicated excess volume.

- 2 Find the optimal blocking and hybridization temperature for your probe design in [Table 9](#).

Table 9 Blocking and Hybridization temperature based on probe design

| Probe Design | Optimal Blocking/Hybridization Temperature |
|---|--|
| SureSelect XT Human All Exon V8/V8+UTR/V8+NCV SureSelect XT Human All Exon V7 SureSelect XT PreCap Clinical Research Exome V4 | 67.5°C |
| All other XT probe designs | 65°C |

- 3 Preprogram a thermal cycler for pre-hybridization blocking as shown in [Table 10](#). Set the heated lid to 105°C. Prewarm the instrument before samples are loaded in [step 5](#).

Table 10 Thermal cycler program for blocking prior to hybridization

| Segment | Number of Cycles | Temperature | Time |
|---------|------------------|---|---------------------------|
| 1 | 1 | 95°C | 5 minutes |
| 2 | 1 | 67.5°C or 65°C (see Table 9) | Hold (at least 5 minutes) |

- To each library sample well, add 5 μL of Blocker Mix, ILM (blue cap). Seal the wells then vortex at high speed for 5 seconds. Spin briefly to collect the liquid and release any bubbles.

CAUTION

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

- Transfer the sample strip to the thermal cycler and run the program in [Table 10](#).

Prepare the additional hybridization reagents described in [step 6](#) and [step 7](#) below during the blocking thermal cycler program. Make sure that the library + Blocker Mix samples are held at 67.5°C or 65°C for at least 5 minutes in Segment 2 before adding the remaining hybridization reaction components in [step 8](#) on [page 15](#).

- Prepare a 25% solution of SureSelect RNase Block according to [Table 11](#). Mix well and keep on ice.

Table 11 Preparation of RNase Block solution (includes excess)

| Reagent | Volume for 1 Hyb | Volume for 8 Hybs (includes excess) | Volume for 24 Hybs (includes excess) |
|-------------------------------------|-------------------|-------------------------------------|--------------------------------------|
| SureSelect RNase Block (purple cap) | 0.5 μL | 4.5 μL | 13.0 μL |
| Nuclease-free water | 1.5 μL | 13.5 μL | 39.0 μL |
| Total | 2 μL | 18 μL | 52 μL |

NOTE

Prepare the mixture described in [step 7](#) below at room temperature at the time of use. Do not keep solutions containing the probe at room temperature for extended periods.

For larger run sizes, hybridization reactions can be started in [step 8](#) below using a multichannel pipette after dividing the master mix volume prepared in [step 7](#) into wells of a strip tube. Using this protocol modification, which includes additional pipetting losses, may require preparation of a larger reagent volume mixture.

- 7 Prepare the Probe Hybridization Mix appropriate for your probe design size. Use [Table 12](#) for probes ≥ 3 Mb or [Table 13](#) for probes < 3 Mb. For custom probes, see the probe tube label for design size range.

Combine the listed reagents at room temperature. Mix well by vortexing at high speed for 5 seconds then spin down briefly. Keep at room temperature until added to wells in [step 8](#).

Table 12 Preparation of Probe Hybridization Mix for **probes ≥ 3 Mb**

| Reagent | Volume for 1 Hyb | Volume for 8 Hybs (includes excess) | Volume for 24 Hybs (includes excess) |
|--|-----------------------------|-------------------------------------|--------------------------------------|
| 25% RNase Block solution (from step 6) | 2 μ L | 18 μ L | 50 μ L |
| Probe (with design ≥ 3 Mb) | 5 μ L | 45 μ L | 125 μ L |
| SureSelect Hybridization Buffer (from step 1) | 13 μ L | 117 μ L | 325 μ L |
| Total | 20 μL | 180 μL | 500 μL |

Table 13 Preparation of Probe Hybridization Mix for **probes < 3 Mb**

| Reagent | Volume for 1 Hyb | Volume for 8 Hybs (includes excess) | Volume for 24 Hybs (includes excess) |
|--|-----------------------------|-------------------------------------|--------------------------------------|
| 25% RNase Block solution (from step 6) | 2 μ L | 18 μ L | 50 μ L |
| Probe (with design < 3 Mb) | 2 μ L | 18 μ L | 50 μ L |
| SureSelect Hybridization Buffer (from step 1) | 13 μ L | 117 μ L | 325 μ L |
| Nuclease-free water | 3 μ L | 27 μ L | 75 μ L |
| Total | 20 μL | 180 μL | 500 μL |

- 8 Keeping the library + Blocker samples in the cycler held at 67.5°C or 65°C, transfer 20 μ L of the room-temperature probe hybridization mix from [step 7](#) to each sample well.
Mix well by pipetting up and down slowly 8 to 10 times.
The hybridization reaction wells now contain approximately 29 μ L.
- 9 Seal the wells with fresh strip caps. Make sure that all wells are completely sealed. Vortex briefly, then spin briefly to remove any bubbles. Immediately return the hybridization strip to the thermal cycler.
- 10 Incubate the hybridization reactions in the thermal cycler for 16 to 24 hours at 67.5°C or 65°C (see [Table 9](#)) with heated lid set to 105°C.

Step 3. Prepare streptavidin beads and buffers for capture

The capture workflow segment uses the components listed in [Table 14](#).

Table 14 Reagents for Capture

| Storage Location | Kit Component | Preparative Steps | Where Used |
|---|--|--|-------------------------|
| +4°C | SureSelect Streptavidin Beads (bottle) | Remove from 4°C just before use, vortex to mix | page 16 |
| SureSelect Max Target Enrichment Kit Overnight Hyb Module Box 1, stored at RT | SureSelect Binding Buffer (bottle) | Ready to use | page 16 |
| | SureSelect Wash Buffer 1 (bottle) | Ready to use | page 17 |
| | SureSelect Wash Buffer 2 (bottle) | Ready to use | page 16 |

NOTE

Set the thermal cycler heated lid to 105°C for all thermal cycler incubation steps used for bead preparation and library capture, including the 6X capture wash incubations at 70°C on [page 17](#).

- 1 Prepare 200 µL aliquots of Wash Buffer 2 for later use in capture wash steps. Using fresh strip tubes, aliquot 6 wells of buffer for each sample in the run. Keep the aliquots at room temperature until directed to pre-warm for washes on [page 17](#).
- 2 Vigorously resuspend the vial of SureSelect Streptavidin Beads on a vortex mixer. The magnetic beads settle during storage.
- 3 For each hybridization sample, add 50 µL of the resuspended beads to wells of a fresh PCR strip tube.
- 4 Equilibrate the beads in SureSelect Binding Buffer using 3× washes:
 - a Add 200 µL of SureSelect Binding Buffer.
 - b Mix by pipetting up and down 20 times or cap the wells and vortex at high speed for 5–10 seconds then spin down briefly.
 - c Put the strip tube into a magnetic separator device.
 - d Wait 5 minutes or until the solution is clear, then remove and discard the supernatant.
 - e Repeat [step a](#) through [step d](#) two more times for a total of 3 washes.
- 5 Resuspend the washed streptavidin beads in 200 µL of SureSelect Binding Buffer.
- 6 Pre-warm the streptavidin capture beads and Binding Buffer by incubating the strip tube in a thermal cycler at 68°C for 10 minutes.

Step 4. Capture the hybridized libraries

- 1 After the 16 to 24 hour hybridization incubation period is complete and after finishing all streptavidin bead preparation steps, add the hybridization mixture to the pre-warmed capture beads as described below.
 - a Using a multichannel pipette, transfer the entire volume (approximately 29 µL) of the hybridization mixtures to wells of pre-warmed beads, while keeping the bead strip in the thermal cycler at 68°C.

- b Using a P200 pipette set to 150 μL , mix by pipetting up and down 8–10 times; make sure the beads are fully resuspended. Seal the wells with fresh caps.
- 2 Incubate the capture reaction strip in the thermal cycler at 68°C for 10 minutes.
- 3 Transfer the SureSelect Wash Buffer 2 aliquots prepared on [page 16](#) (200 μL /well) to a thermal cycler held at 70°C. Pre-warm the buffer until used for washes in [step 8](#).
- 4 When the 10 minute capture incubation ([step 2](#)) is complete, spin the samples briefly to collect the liquid.
- 5 Put the capture reaction strip in a magnetic separator to collect the beads. Wait until the solution is clear (approximately 1 to 2 minutes), then remove and discard the supernatant.
- 6 Resuspend the beads in 200 μL of SureSelect Wash Buffer 1. Mix by pipetting up and down 15–20 times, until beads are fully resuspended.
- 7 Return the sample strip to the magnetic separator. Wait for the solution to clear (approximately 1 minute), then remove and discard the supernatant.

Spin the samples briefly to collect any residual wash buffer and remove with a P20 pipette.

CAUTION

To ensure specificity of capture, make sure that the SureSelect Wash Buffer 2 is pre-warmed to 70°C before use in the wash steps below.

During the washes in [step 8](#), keep the sample strip at room temperature during the pipetting and vortexing steps prior to each 70°C incubation step.

Do not use a tissue incubator, or other devices with significant temperature fluctuations, for the 70°C incubation steps.

-
- 8 **Place the capture sample strip in a rack at room temperature.** Wash the beads with Wash Buffer 2, using the steps below.
 - a Resuspend the beads in 200 μL of 70°C pre-warmed Wash Buffer 2. **Make sure the beads are completely in suspension before proceeding.** To ensure thorough resuspension, Agilent recommends the following two-part mixing routine. Keep the capture reactions at room temperature while mixing.
 - i) Using a P200 pipette set to 150 μL , pipette up and down 15–20 times, until beads are fully resuspended. Pipette slowly and gently to reduce bubble formation.
 - ii) Seal the wells with fresh caps and vortex at high speed for 8 seconds. Spin the strip tube briefly to collect the liquid and release any bubbles without pelleting the beads.
 - b Incubate the samples for 5 minutes at 70°C in a thermal cycler.
 - c Put the sample strip in the magnetic separator at room temperature.
 - d Wait 1 minute for the solution to clear, then remove and discard the supernatant. Return samples to the rack at room temperature.
 - e Repeat [step a](#) through [step d](#) five more times for a total of 6 washes.

When the final wash is complete, spin the samples briefly to collect any residual Wash Buffer 2 and remove with a P20 pipette.

- 9 Add 24 μL of nuclease-free water to each sample well. Pipette up and down 8 times to resuspend the beads.
- 10 Keep the samples on ice until they are used in the PCR reactions on [page 19](#).

NOTE

The captured DNA or cDNA libraries are retained on the streptavidin beads during the post-capture amplification step.

Step 5. Amplify the captured libraries

The post-capture amplification workflow segment uses the components listed in [Table 15](#). Prepare as directed before use (refer to the *Where Used* column).

Table 15 Reagents for post-capture amplification

| Storage Location | Kit Component | Preparative Steps | Where Used |
|--|--|--|-------------------------|
| SureSelect Max Target Enrichment Kit Overnight Hyb Module Box 2, stored at -20°C | Amplification Master Mix (red cap or bottle) | Thaw on ice then keep on ice. Mix thoroughly by inversion at least 5X. Do not vortex. | page 19 |
| SureSelect Max Blockers and Primers Module for ILM, stored at -20°C | SureSelect Post-Capture Primer Mix (clear cap) | Thaw and keep on ice, vortex to mix. | page 19 |
| +4°C | SureSelect Max Purification Beads OR AMPure XP Beads | Equilibrate at room temperature for at least 30 minutes before use, vortex to mix. | page 20 |

- 1 Preprogram a thermal cycler as shown in [Table 16](#). Set the heated lid to 105°C. Prewarm the instrument before samples are loaded in [step 6](#) on [page 19](#).

Table 16 Post-Capture PCR thermal cycler program (50 µL vol)

| Segment | Number of Cycles | Temperature | Time |
|---------|--|-------------|------------|
| 1 | 1 | 98°C | 45 seconds |
| 2 | 10–16 (see Table 17 for probe design size-based cycle number recommendation) | 98°C | 15 seconds |
| | | 60°C | 30 seconds |
| | | 72°C | 30 seconds |
| 3 | 1 | 72°C | 1 minute |
| 4 | 1 | 4°C | Hold |

Table 17 Post-capture PCR cycle number recommendations

| Probe Design Size | Cycles | |
|--|---------------------|---------------------|
| | DNA input libraries | RNA input libraries |
| Probes <0.2 Mb | 16 cycles | 16 cycles |
| Probes 0.2–3 Mb | 12–16 cycles | 14 cycles |
| Probes 3–5 Mb | 11–12 cycles | 13 cycles |
| Probes >5 Mb (including Human All Exon and Exome probes) | 10–11 cycles | 12 cycles |

NOTE

To avoid cross-contamination, it is recommended to set up PCR reactions (all components except the libraries) in a dedicated clean area or PCR hood with UV sterilization and positive air flow.

- 2 Mix the thawed Amplification Master Mix (red cap or bottle) thoroughly by inversion (do not vortex) then spin briefly.
- 3 Prepare the appropriate volume of post-capture PCR reaction mix as described in [Table 18](#), on ice. Mix well on a vortex mixer.

Table 18 Preparation of post-capture PCR reaction mix

| Reagent | Volume for 1 reaction | Volume for 8 reactions (includes excess) | Volume for 24 reactions (includes excess) |
|--|-----------------------------|--|---|
| Amplification Master Mix (red cap or bottle) | 25 μ L | 225 μ L | 650 μ L |
| SureSelect Post-Capture Primer Mix (clear cap) | 1 μ L | 9 μ L | 26 μ L |
| Total | 26 μL | 234 μL | 676 μL |

- 4 Add 26 μ L of the PCR reaction mix prepared in [Table 18](#) to each sample well containing 24 μ L of bead-bound target-enriched library DNA or cDNA.
- 5 Mix the PCR reactions well by pipetting up and down until the bead suspension is homogeneous. Avoid splashing samples onto well walls; do not spin the samples at this step.
- 6 Place the sample strip in the thermal cycler and run the program in [Table 16](#).

NOTE

The SureSelect Max Purification Beads or AMPure XP Beads used in the next section must be equilibrated to room temperature for at least 30 minutes before use.

- 7 When the PCR amplification program is complete, spin the amplified library strip briefly. Proceed to purification.

NOTE

In the following purification protocol, the hybrid capture streptavidin beads remain mixed with the magnetic purification beads throughout the procedure.

If preferred, you can remove the streptavidin beads prior to purification by adding the following steps: 1) Place the amplified library strip tube on the magnetic stand at room temperature 2) Wait 2 minutes for the solution to clear 3) Transfer each supernatant (approximately 50 μ L) to fresh strip tube wells and proceed with the purification procedure. The streptavidin beads can be discarded after collecting the library solution.

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 the final libraries using magnetic purification beads

Purify the amplified libraries using room-temperature SureSelect Max Purification Beads or AmpPure XP Beads. Critical purification protocol parameters are summarized for experienced users in [Table 19](#).

Table 19 Magnetic purification bead cleanup parameters after post-capture PCR

| Parameter | Value |
|---|--------------------------|
| Volume of RT purification bead suspension added to each sample well | 50 μ L |
| Final elution solvent and volume | 25 μ L Low TE Buffer |
| Amount of eluted sample transferred to fresh well | Approximately 24 μ L |

- 1 Prepare 400 μ 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 Add 50 μ L of the purification bead suspension to each amplified library sample well, containing approximately 50 μ L of streptavidin bead suspension.
- 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 strip tube into a magnetic separation device. Wait for the solution to clear (approximately 2 to 5 minutes).
- 7 Keep the 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 strip tube in the magnetic stand while you dispense 200 μ 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 strip tube at room temperature for 2–5 minutes or until the residual ethanol has just evaporated.
- 13 Elute the library by adding 25 μ L of 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 or cDNA fragments.
- 16 Put the strip tube in the magnetic stand and leave until the solution is clear (up to 5 minutes).

17 Transfer the cleared supernatant (approximately 24 µL) to a fresh well. Keep on ice.

You can discard the beads at this time.

NOTE

If magnetic beads are carried over during the final elution step, samples can be placed on the magnetic stand while removing an aliquot for QC (below) and while removing an aliquot for pooling on [page 23](#).

Stopping Point If you do not plan to continue through the library pooling for NGS step on same day, seal the wells and store at 4°C overnight or at –20°C for prolonged storage (remove aliquot for QC analysis before storage, if appropriate).

Step 7. QC and quantify final libraries

Analyze a sample of each library using one of the platforms listed in [Table 20](#). Follow the instructions in the linked user guide provided for each assay.

Table 20 Post-capture 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 | High Sensitivity D1000 ScreenTape | Agilent High Sensitivity D1000 Assay Quick Guide | 2 µL |
| Agilent 5200/5300/5400 Fragment Analyzer system | HS NGS Fragment Kit (1-6000 bp) | Agilent HS NGS Fragment Kit (1-6000 bp) Kit Guide | 2 µL |

Each analysis method provides an electropherogram showing the size distribution of fragments in the sample and tools for determining the concentration of library DNA or cDNA in the sample. 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 concentration. See [Table 21](#) for fragment size distribution guidelines.

Table 21 Post-capture expected library fragment size guidelines

| Hybridization input | Expected average fragment size (150–1000 bp region) |
|--|---|
| DNA library prepared from intact DNA fragmented for 2x100 NGS | 350 to 450 bp |
| DNA library prepared from intact DNA fragmented for 2x150 NGS | 380 to 480 bp |
| DNA library prepared from FFPE DNA fragmented for 2x100 or 2x150 NGS | 250 to 390 bp |
| cDNA library prepared from intact RNA | 380 to 480 bp |
| cDNA library prepared from FFPE RNA | 250 to 390 bp |

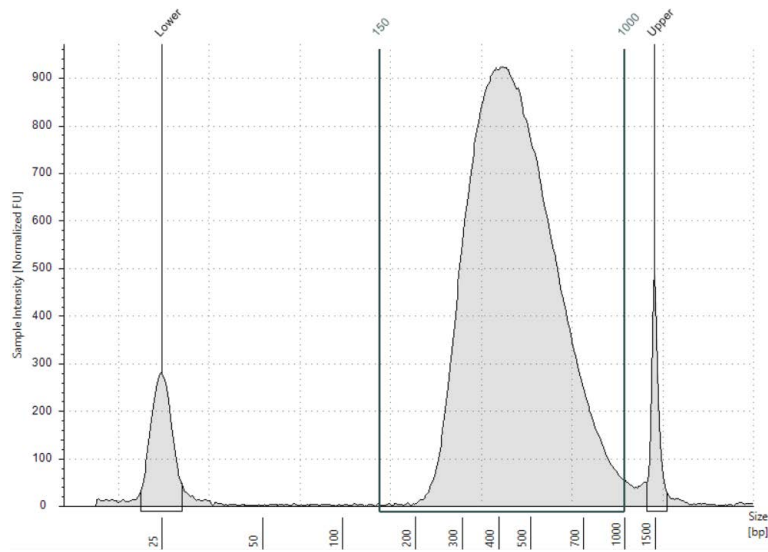


Figure 2 Post-capture library prepared from a high-quality gDNA sample analyzed using a High Sensitivity D1000 ScreenTape assay.

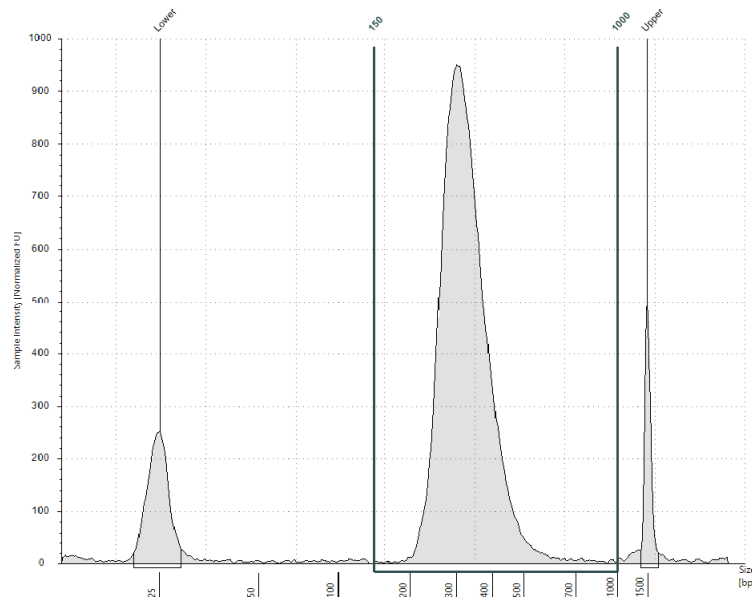


Figure 3 Post-capture library prepared from an FFPE gDNA sample analyzed using a High Sensitivity D1000 ScreenTape assay.

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

Post-capture pooling guidelines

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.

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.

Combine the library samples or pre-capture pool 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 or cDNA 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 22 shows an example of the amount of 4 index-tagged samples (of different concentrations) and Low TE needed for a final volume of 20 μL at 10 nM DNA or cDNA.

Table 22 Example of volume calculation for total volume of 20 μL at 10 nM concentration

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

Store the library pool under the conditions specified by your sequencing provider. Typically, libraries should be stored at -20°C for short-term storage.

3

Appendix: NGS and Analysis Guidelines

| | |
|---|----|
| SureSelect Max Library Composition | 25 |
| Sequencing Setup and Run Guidelines | 26 |
| Analysis Pipeline Guidelines | 27 |
| AGeNT software guidelines for SureSelect Max DNA or RNA workflows | 28 |
| RNA strandedness guidelines | 28 |
| SureSelect Max UDI Information | 28 |
| SureSelect Max index sequences | 30 |

This appendix provides guidelines for completing NGS using the Illumina platform and for SureSelect Max library read processing steps.

SureSelect Max Library Composition

The final SureSelect Max library fragments prepared using Agilent's Illumina-compatible kit modules are shown in [Figure 4](#). 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.

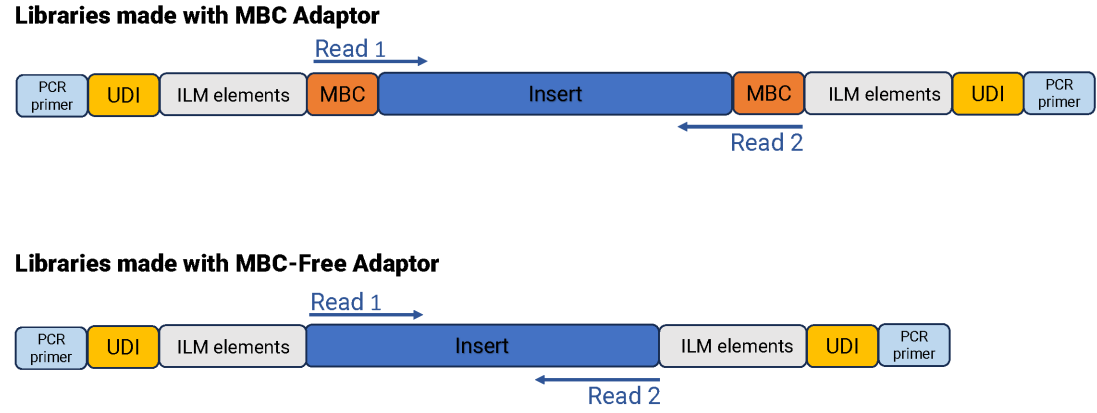


Figure 4 Content of SureSelect Max for ILM sequencing libraries. Each fragment contains one target insert (dark blue) surrounded by the following elements: Illumina paired-end sequencing elements (gray), unique dual indexes (gold), library PCR primers (light blue) and optional molecular barcode (MBC) sequences. MBCs include a 3-bp degenerate barcode plus dark bases.

Sequencing Setup and Run Guidelines

Proceed to cluster amplification using the appropriate Illumina Paired-End Cluster Generation Kit. [Table 23](#) provides guidelines for use of several Illumina instrument and chemistry combinations suitable for processing the SureSelect Max target-enriched NGS libraries. For other Illumina NGS platforms, consult Illumina’s documentation for kit configuration and seeding concentration guidelines.

Table 23 Illumina Kit Configuration Selection Guidelines

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

Seeding concentration and cluster density may also need to be optimized based on the fragment size range for the library and on the desired output and data quality. Begin optimization using a seeding concentration in the middle of the range listed in [Table 23](#) 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 24](#) showing example settings for 2x150 bp sequencing.

Table 24 Run settings for 2x150 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 28](#).
- 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 MBC-tagged libraries, turn off any adaptor trimming tools included in Illumina's run setup and read processing software applications. Adaptors are trimmed in later processing steps to ensure proper processing of the degenerate MBC sequences in the adaptors.
- 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. For use in these applications, the SureSelect Max index sequences can be obtained by downloading the [SureSelect Max Index Sequence Resource](#) Excel spreadsheet from Agilent.com. The provided sequences should be converted to .tsv/.csv file format or copied to a Sample Sheet according to Illumina's specifications for each application. If you need assistance with SureSelect Max run setup in your selected application (e.g., generating index files or Sample Sheet templates), 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 appropriate for SureSelect Max DNA and RNA libraries. Your NGS pipeline may vary.

- Demultiplex using Illumina's bcl2fastq, BCL Convert or DRAGEN software to generate paired end reads based on the dual indexes and remove sequences with incorrectly paired P5 and P7 indexes. For MBC-tagged libraries, 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. For MBC-free libraries, trimming tools in Illumina's demultiplexing software can be turned on.
- If you have prepared MBC-tagged libraries, but your sequence analysis pipeline excludes MBCs, you can remove MBCs during the demultiplexing step by trimming or masking the first five bases from each read. See [page 41](#) for guidelines.
- For MBC-tagged reads, the demultiplexed FASTQ data needs to be pre-processed to trim the sequencing adaptors and to extract and use MBC sequences for de-duplication. Agilent's AGeNT toolkit, described on [page 28](#), can be used for these pre-processing steps.

NOTE

Read pre-processing steps can also be completed using suitable open-source software tools, such as fgbio. Performance of open-source tools should be verified for appropriate adaptor and MBC sequence processing on both strands. Some non-Agilent adaptor trimmers may fail to remove the MBC sequences from the opposite adaptor, which may affect alignment quality.

AGeNT software guidelines for SureSelect Max DNA or RNA workflows

Agilent's AGeNT is a Java-based toolkit for library read processing steps, designed for users with bioinformatics expertise to enable building internal analysis pipelines. To download this toolkit, visit the [AGeNT page at www.agilent.com](http://www.agilent.com). Use of the AGeNT read processing tools is outlined briefly below. See the [AGeNT Best Practices](#) document for more information. Instructions provided for the SureSelect XT HS2 platform are also applicable to the SureSelect Max platform.

- Prior to variant discovery, demultiplexed SureSelect Max library FASTQ data are pre-processed to remove sequencing adaptors and extract the MBC sequences (if present) using the AGeNT Trimmer module.
- The trimmed reads should be aligned (and MBC tags added to the aligned BAM file where applicable) using a suitable tool such as BWA-MEM for DNA libraries or STAR for RNA libraries.
- For MBC-containing libraries only, the AGeNT CReaK (Consensus Read Kit) tool is used to generate consensus reads and mark or remove duplicates for the aligned reads based on the read MBCs. CReaK includes consensus generation options suitable for DNA libraries (duplex, hybrid, or single-strand mode) and RNA libraries (single-strand mode).

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

RNA strandedness guidelines

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

SureSelect Max UDI Information

The SureSelect Max unique dual indexes (UDIs) are added to the library fragments during the library preparation workflow segment. Each fragment contains a unique 8-bp P5 and P7 index (see [Figure 4](#) on page 25) suitable for Illumina sequencing platforms.

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.

Index sequences are provided on [page 30](#) through [page 37](#). 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 26](#) through [Table 33](#) 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 two orientations (forward and reverse complement) for use with different platforms and sequencing run setup and management tools, e.g., Local Run Manager and Instrument Run Setup. A selection of Illumina sequencing platforms and their P5 sequencing orientations are shown in [Table 25](#). Correct representation of the P5 index orientation in sample sheets or during sequencing run setup is crucial to successful demultiplexing. Refer to Illumina support documentation and resources to determine the correct P5 index orientation for your application.

Table 25 P5 index sequencing orientation by Illumina platform

| P5 Index Orientation | Platform |
|----------------------|--|
| Forward | MiSeq |
| Reverse Complement* | NovaSeq 6000 with v1.5 chemistry NextSeq 500/550/1000/2000 iSeq 100 MiniSeq |

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

SureSelect Max index sequences

Index sequences can also be obtained by downloading the [SureSelect Max Index Sequence Resource](#) Excel spreadsheet from Agilent.com.

Table 26 SureSelect Max Index Primer Pairs 1–48, provided in orange 96-well plate or in strip tubes

| Primer Pair # | Index Strip | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement | Primer Pair # | Index Strip | 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 | GTTCTCAT |
| 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 | GTA ACTTG |
| 24 | H03 | TGGCACCA | TGCTGCTC | GAGCAGCA | 48 | H06 | CGCGCAAT | CAGAGCAG | CTGCTCTG |

Table 27 SureSelect Max Index Primer Pairs 49–96, provided in orange 96-well plate

| Primer Pair # | Index Strip | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement | Primer Pair # | Index Strip | 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 28 SureSelect Max Index Primer Pairs 97–144, provided in blue 96-well plate

| Primer Pair # | Index Strip | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement | Primer Pair # | Index Strip | 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 | TGCGATTC |
| 120 | H03 | AGACGCCT | GCGCAATT | AATTGCGC | 144 | H06 | CTCCGTT | CTTCACGT | ACGTGAAG |

Table 29 SureSelect Max Index Primer Pairs 145–192, provided in blue 96-well plate

| Primer Pair # | Index Strip | P7 Index Forward | P5 Index Forward | P5 Index Reverse Complement | Primer Pair # | Index Strip | 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 | GTAATTCA |
| 152 | H07 | CACATTCG | TGCAATGC | GCATTGCA | 176 | H10 | TGATTGGC | CGAATCAT | ATGATTCCG |
| 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 | TGTTCCGG |
| 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 30 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 | GTTAGTT | 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 | GACTGATG | 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 31 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 32 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 | AGGTCAT | CGAGTCGA | TCGACTCG |
| 303 | G02 | CGGTCTCT | GAGATCGC | GCGATCTC | 327 | G05 | GACGTGCA | AGGTCAT | 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 33 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 |

4 Reference

Reagent Kit Contents [39](#)

Troubleshooting Guide [40](#)

Quick Reference Protocol: Max Overnight Hyb Target Enrichment [42](#)

This section contains reference information, including Reagent Kit contents, troubleshooting information and a quick-reference protocol for experienced users.

Reagent Kit Contents

SureSelect Max Target Enrichment with overnight hybridization workflow uses the kits listed in [Table 34](#). Detailed contents of each of the multi-part component kits are shown in [Table 35](#) through [Table 37](#).

Table 34 Kits for SureSelect Max Target Enrichment with Overnight Hybridization

| SureSelect Max Kits and Included Component Kits | Component Kit Part Numbers | | Storage Condition |
|---|----------------------------|-------------------|-------------------|
| | 16 Hybridizations | 96 Hybridizations | |
| SureSelect Max Overnight Hyb Kit (G9690A/G9690B) | | | |
| SureSelect Max Target Enrichment Kit Overnight Hyb Module, Box 1 | 5282-0132 | 5282-0134 | Room Temperature |
| SureSelect Max Target Enrichment Kit Overnight Hyb Module, Box 2 | 5282-0133 | 5282-0135 | -20°C |
| SureSelect Streptavidin Beads | 5191-5741 | 5191-5742 | +4°C |
| SureSelect Max Blockers and Primers Module for ILM (G9699A/G9699B) | 5282-0136 | 5282-0137 | -20°C |
| SureSelect Max Purification Beads (G9962A/G9962B) | 5282-0225 | 5282-0226 | +4°C |

Component Kit Details

Table 35 SureSelect Max Target Enrichment Kit Overnight Hyb Module, Box 1 content

| Kit Component | 16 Hyb Kit (p/n 5282-0132) | 96 Hyb Kit (p/n 5282-0134) |
|---------------------------|----------------------------|----------------------------|
| SureSelect Hyb 1 | tube with orange cap | bottle |
| SureSelect Hyb 2 | tube with red cap | tube with red cap |
| SureSelect Hyb 4 | tube with black cap | tube with black cap |
| SureSelect Binding Buffer | bottle | bottle |
| SureSelect Wash Buffer 1 | bottle | bottle |
| SureSelect Wash Buffer 2 | bottle | bottle |

Table 36 SureSelect Max Target Enrichment Kit Overnight Hyb Module, Box 2 content

| Kit Component | 16 Hyb Kit (p/n 5282-0133) | 96 Hyb Kit (p/n 5282-0135) |
|--------------------------|----------------------------|----------------------------|
| SureSelect Hyb 3 | tube with yellow cap | tube with yellow cap |
| SureSelect RNase Block | tube with purple cap | tube with purple cap |
| Amplification Master Mix | tube with red cap | bottle |

Table 37 SureSelect Max Blockers and Primers Module for ILM content

| Kit Component | 16 Hyb Kit (p/n 5282-0136) | 96 Hyb Kit (p/n 5282-0137) |
|------------------------------------|----------------------------|----------------------------|
| Blocker Mix, ILM | tube with blue cap | tube with blue cap |
| SureSelect Post-Capture Primer Mix | tube with clear cap | tube with clear cap |

Troubleshooting Guide

If vacuum concentrator is not available for library volume reduction prior to hybridization

- ✓ The standard hybridization protocol requires addition of prepared library samples or library pools in a 4 μL volume, using a vacuum concentrator to reduce the volume of the library pools where needed. If a vacuum concentrator is not available in your laboratory, use the protocol modifications below to generate concentrated library pools for hybridization. This modification may result in some loss of library complexity.

Post-capture pooling workflow:

- Method 1: During the SureSelect Max Library Preparation protocol using Enzymatic Fragmentation or Covaris shearing, in the final pre-capture library preparation clean-up step reduce the amount of nuclease-free water used for elution from 15 μL to 6 μL .
- Method 2: If prepared libraries were already eluted in 15 μL , libraries can be concentrated using an additional round of magnetic bead purification. Follow the purification instructions on [page 20](#), modifying [step 3](#) to use a bead volume of 1.8X the available sample volume, then elute using 6 μL of nuclease-free water in [step 13](#).

Pre-capture pooling workflow:

- For hybridization strip setup on [page 12](#), make pools containing 2.5X amount of each library given in [Table 7](#) on page 12. For example, for gDNA libraries enriched with the SureSelect XT PreCap Human All Exon V8 probe, pool 1250 ng of each of 8 indexed libraries for total pool of 10 μg library DNA.
- Concentrate each pool using an additional round of magnetic bead purification. Follow the purification instructions on [page 20](#), modifying [step 3](#) to use a bead volume of 1.8X the total sample volume.
- Elute in 8 μL of nuclease-free water in [step 13](#). Place 4 μL of each concentrated pool in a separate well for the hybridization reaction. The remaining concentrated pool volume is not used in the hybridization reaction.

If yield of captured libraries is low

- ✓ The protocol includes specific thawing, temperature control, pipetting, and mixing instructions which are required for optimal performance. Be sure to adhere to all instructions when setting up the reactions.
- ✓ PCR cycle number may require optimization. Repeat target enrichment for the sample, increasing the post-capture PCR cycle number by 1 to 2 cycles.
- ✓ The probe used for hybridization may have been compromised. Verify the expiration date on the probe vial or Certificate of Analysis. Adhere to the recommended storage and handling conditions. Ensure that the probe hybridization mix is prepared at the time of use, as directed on [page 15](#), and that solutions containing the probe are not held at room temperature for extended periods.
- ✓ 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](#)). 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 or cDNA fragments.

If captured library fragment size is different than expected in electropherograms

- ✓ Libraries prepared from FFPE DNA or RNA samples may have a smaller fragment size distribution due to the presence of fragments in the sample input that are smaller than the optimal fragment size for target enrichment. Adhere to the FFPE DNA or RNA quality guidelines provided in the SureSelect Max Library Preparation Module user guides.
- ✓ Library fragment size selection during post-capture 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 and verify that you are using the bead volume recommended for post-capture purification on [page 20](#).

If low percent on target is observed in library sequencing results

- ✓ Stringency of post-hybridization washes may have been lower than required. Complete the wash steps as directed, paying special attention to the details of SureSelect Wash Buffer 2 washes listed below:
 - Ensure that SureSelect Wash Buffer 2 is pre-warmed to 70°C before use (see [page 17](#)). Select a thermal cycler with a block configured for efficient heating of 0.2 mL liquid volumes; ensure that the plasticware containing the wash buffer is fully seated in the thermal cycler block wells, with minimal liquid volume visible above the block during the pre-warming step.
 - Bead suspensions are mixed thoroughly during washes by pipetting up and down and vortexing (see [page 17](#)).
- ✓ Minimize the amount of time that hybridization reactions are exposed to RT conditions during hybridization setup. Locate a vortex and plate spinner or centrifuge in close proximity to thermal cycler to retain the elevated sample temperature during mixing and transfer steps ([step 8](#) to [step 10](#) on [page 15](#)).

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 using one of the options below:
 - If demultiplexing using bcl2fastq, MBCs may be masked by including the base mask **N5Y*,I8,I8,N5Y***. If demultiplexing using BCL Convert, MBCs may be trimmed by including the following string in the sample sheet header: **OverrideCycles,N5Y*,I8,I8,N5Y***. For both methods, * is replaced with value equal to the remaining read length after masking or trimming. For example, use **N5Y146,I8,I8,N5Y146** for 2x150 NGS with 151 cycles (as shown in [Table 24](#) on page 26). The sum of the values following N and Y must match the read length value in the RunInfo.xml file.
 - The first 5 bases may be trimmed from the demultiplexed FASTQ files using the AGeNT Trimmer module while trimming adaptor sequences or using a suitable processing tool of your choice, such as seqtk. Non-Agilent adaptor trimmer performance should be verified for removal of the MBC sequences from the opposite adaptor (refer to [Figure 4](#) on page 25); failure to remove MBC sequences from both strands may affect alignment quality.

Quick Reference Protocol: Max Overnight Hyb Target Enrichment

An abbreviated protocol summary is provided below for experienced users. Use the complete protocol on [page 10](#) to [page 21](#) until you are familiar with all of the protocol details.

| Step | Summary of Conditions |
|--|---|
| Hybridization (Day 1) | |
| Place libraries in hyb wells and adjust volumes to 4 µL/well | Place prepared libraries or library pools into strip tube wells and adjust volume to 4 µL as below. For post-cap pooling: Use 12 µL of undiluted prepared library (recommended) or use at least 500 ng gDNA library or 200 ng cDNA library per well. For pre-cap pooling: Pool libraries into wells according to Table 7 on page 12. Reduce well volumes to <4 µL by vacuum concentration at ≤45°C. Bring well volumes to 4 µL with nuclease-free H ₂ O> vortex> spin> keep on ice. |
| Prep Overnight Hyb Buffer | Per 8 Hyb reactions: 66.3 µL Hyb 1 + 2.7 µL Hyb 2 + 26.5 µL Hyb 3 + 34.5 µL Hyb 4 Per 24 reactions: 185.6 µL Hyb 1 + 7.6 µL Hyb 2 + 74.2 µL Hyb 3 + 96.6 µL Hyb 4 Prepare at RT> mix> spin> keep at RT. |
| Program thermal cycler | Input the Blocking/Hybridization program in Table 38 , using the temperature appropriate for your probe in segment 2. Pre-warm before loading samples. |
| Run pre-hybridization blocking protocol | Add 5 µL Blocker Mix ILM to each sample well; vortex> spin> place strip in cycler> resume program. Block at 67.5°C or 65°C for at least 5 min while preparing the remaining hyb reagents. |
| Prepare 25% RNase Block dilution | Per 8 Hyb reactions: 4.5 µL RNase Block + 13.5µL nuclease-free H ₂ O Per 24 reactions: 13 µL RNase Block + 39 µL nuclease-free H ₂ O Prepare on ice> mix> spin> keep on ice. |
| Prepare Probe/Hyb Mix | Prepare the Probe/Hyb Mix for your probe design size--see Table 39 . Prepare at RT> vortex> spin> keep briefly at RT. |
| Run the hybridization | Keep samples in cycler at 67.5°C/65°C while adding 20 µL Probe/Hyb Mix to wells> mix by pipetting. Seal wells completely> vortex> spin> return strip to cycler> hybridize 16 to 24 hours. |

Table 38 Blocking/Hybridization program (29 µL vol; heated lid at 105°C)

| Segment | Probe | Number of Cycles | Temperature (Purpose) | Time |
|---------|--|------------------|-----------------------|---|
| 1 | All probes | 1 | 95°C (Pre-blocking) | 5 minutes |
| 2 | SureSelect XT Human All Exon V8/V8+UTR/V8+NCV, Human All Exon V7 or Clinical Research Exome V4 | 1 | 67.5°C (Blocking/Hyb) | Hold: Do blocking for ≥5 min Do hybridization for 16–24 hours |
| | All other SureSelect XT probe designs | 1 | 65°C (Blocking/Hyb) | Hold: Do blocking for ≥5 min Do hybridization for 16–24 hours |

Table 39 Preparation of Probe/Hyb Mix

| Reagent | Probe designs ≥3 Mb | | | Probe designs <3 Mb | | |
|-------------------------------|---------------------|---------------|---------------|---------------------|---------------|---------------|
| | Per Hyb Reaction | 8 Hybs | 24 Hybs | Per Hyb Reaction | 8 Hybs | 24 Hybs |
| 25% RNase Block solution | 2 µL | 18 µL | 50 µL | 2 µL | 18 µL | 50 µL |
| Probe | 5 µL | 45 µL | 125 µL | 2 µL | 18 µL | 50 µL |
| Prepared Overnight Hyb Buffer | 13 µL | 117 µL | 325 µL | 13 µL | 117 µL | 325 µL |
| Nuclease-free water | — | — | — | 3 µL | 27 µL | 75 µL |
| Total | 20 µL | 180 µL | 500 µL | 20 µL | 180 µL | 500 µL |

| Step | Summary of Conditions |
|---|--|
| Capture (Day 2) | |
| Aliquot Wash Buffer 2 | Aliquot 6 × 200 µL of SureSelect Wash Buffer 2 for each sample |
| Prepare streptavidin beads | For each sample, wash 50 µL SureSelect Streptavidin beads 3× in 200 µL SureSelect Binding Buffer. After final wash, resuspend beads in 200 µL SureSelect Binding Buffer and pre-warm beads/buffer at 68°C for 10 minutes. |
| Capture hybridized libraries | Add hybridized samples (~29 µL) to pre-warmed beads (200 µL) held in cycler> mix well by pipetting. Incubate capture strip in thermal cycler at 68°C for 10 minutes. During incubation, pre-warm the 6 × 200 µL aliquots of SureSelect Wash Buffer 2 to 70°C. |
| Wash captured libraries | Spin capture strip briefly> collect beads using a magnetic stand> discard supernatant. Wash beads 1× with 200 µL SureSelect Wash Buffer 1 at RT. Wash beads 6× with 200 µL pre-warmed SureSelect Wash Buffer 2 (for each wash: add 200 µL 70°C Wash Buffer 2> mix well by pipetting and vortexing at RT> incubate 5 minutes at 70°C> collect beads at RT> discard supernatant). Resuspend washed beads in 24 µL nuclease-free H ₂ O> pipette to mix> keep on ice. |
| Post-capture amplification (Day 2) | |
| Program thermal cycler | Input the appropriate thermal cycler program in Table 40 . Pre-warm before loading samples. |
| Prepare Post-capture PCR Mix | Per 8 Hyb reactions: 225 µL Amplification Master Mix + 9 µL SureSelect Post-Capture Primer Mix Per 24 reactions: 650 µL Amplification Master Mix + 26 µL SureSelect Post-Capture Primer Mix Prepare on ice> mix> spin> keep on ice. |
| Amplify the bead-bound captured libraries | Add 26 µL Post-capture PCR Mix to each well containing 24 µL captured library bead suspension; pipette to mix (do not spin)> place strip in cycler> run thermal cycler program. |
| Purify amplified libraries | 50 µL amplified library bead suspension + 50 µL purification bead suspension> mix> incubate 5 min> collect beads> discard supernatant. Wash beads 2X with 70% ethanol. Elute in 25 µL Low TE Buffer> mix> incubate 2-5 minutes> collect beads> retain supernatant. |
| Quantify and qualify libraries | Analyze quantity and quality using TapeStation or Fragment Analyzer System |

Table 40 Post-Capture PCR thermal cycler program (50 µL vol; heated lid at 105°C)

| Segment | Number of Cycles | Temperature | Time | |
|---------|--|--|----------------------|--|
| 1 | 1 | 98°C | 45 seconds | |
| 2 | gDNA libraries (DNA input): Probes <0.2 Mb: 16 cycles Probes 0.2–3 Mb: 12–16 cycles Probes 3–5 Mb: 11–12 cycles Probes >5 Mb: 10–11 cycles | OR cDNA libraries (RNA input): Probes <0.2 Mb: 16 cycles Probes 0.2–3 Mb: 14 cycles Probes 3–5 Mb: 13 cycles Probes >5 Mb: 12 cycles | 98°C 60°C 72°C | 15 seconds 30 seconds 30 seconds |
| 3 | 1 | 72°C | 1 minute | |
| 4 | 1 | 4°C | Hold | |

In This Book

This guide provides instructions for SureSelect Max Target Enrichment using an Overnight Hybridization workflow including either post-capture or pre-capture pooling of NGS libraries.

