

HaloPlex Target Enrichment System

For Illumina Sequencing

Protocol

Version H0, February 2025

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



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2015, 2018, 2025

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

G9900-90001

Edition

Version H0, February 2025

Printed in USA

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

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 describes an optimized protocol for using the HaloPlex target enrichment system to prepare sequencing library samples for Illumina paired-end multiplexed sequencing platforms.

1 Before You Begin

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

2 Sample Preparation

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

3 Reference

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

What's New in Version H0

- Support for kits supplied with re-Albumin Solution, replacing BSA Solution (see [page 16](#), [page 17](#), and [page 44](#)). The protocols in this publication are compatible with use of either form of albumin-containing solution.
- Updates to [Technical Support](#) contact information on [page 2](#)
- Updates to [Notices](#) on [page 2](#)
- Updates to required materials supplier information in [Table 1](#) on page 8 and [Table 3](#) on page 10
- Updates to lists of available HaloPlex kits in [Table 2](#) on page 9 and [Table 11](#) on page 43
- Removal of support information for ClearSeq pre-designed probes throughout this publication
- Updates to information on use of Agilent's SureDesign for access to HaloPlex probe designs and kit ordering information on [page 9](#) and [page 12](#)
- Removal of recommendation for hybridization using the SureCycler thermal cycler (see [Table 3](#) on page 10 and [page 27](#))
- Updates to purchasing information for the Agilent 2100 Bioanalyzer instrument in [Table 4](#) on page 11, and related updates to validation instructions on [page 22](#) to [page 23](#) and [page 38](#) to [page 39](#)
- Updates to Agilent.com webpage hyperlinks
- Updates to downstream sequencing support information on [page 41](#)

Content

1 Before You Begin

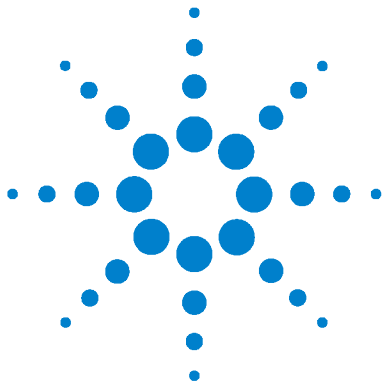
Procedural Notes	7
Safety Notes	7
Required Reagents	8
Required Equipment	10
Optional Validation Reagents and Equipment	11

2 Sample Preparation

Run Size Considerations	14
Run Time Considerations	14
Step 1. Digest genomic DNA with restriction enzymes	15
Step 2. Hybridize digested DNA to HaloPlex probe for target enrichment and sample indexing	25
Step 3. Capture the target DNA	28
Step 4. Ligate the captured, circularized fragments	31
Step 5. Prepare the PCR Master Mix	32
Step 6. Elute captured DNA with NaOH	33
Step 7. PCR amplify the captured target libraries	34
Step 8. Purify the amplified target libraries	35
Step 9. Validate enrichment and quantify enriched target DNA	37
Step 10. Pool indexed samples and perform multiplexed sequencing	41

3 Reference

Kit Contents	43
Nucleotide Sequences of HaloPlex Indexes	46
Qualitative analysis of enrichment by gel electrophoresis	47



1 Before You Begin

Procedural Notes	7
Safety Notes	7
Required Reagents	8
Required Equipment	10
Optional Validation Reagents and Equipment	11

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

Procedural Notes

- The HaloPlex protocol is optimized for digestion of 200 ng of genomic DNA (split among 8 different restriction digestion reactions) plus 25 ng excess DNA, for a total of 225 ng genomic DNA. Using lower amounts of DNA in the enrichment protocol can adversely affect your results. Use a fluorometry-based DNA quantitation method, such as PicoGreen stain or Qubit fluorometry to quantify the DNA starting material.
- Always keep pre-amplification and post-amplification DNA samples in separate work areas. Perform the enrichment procedure in the pre-amplification area. Open and store the amplified, enriched DNA samples only in the post-amplification area.
- Possible stopping points, where DNA samples may be stored between steps, are marked in the protocol. Store the samples at -20°C , but do not subject the samples to multiple freeze/thaw cycles.
- Ensure that master mixes are thoroughly mixed, by pipetting up-and-down or by gentle vortexing, before distributing to the samples.
- In general, follow Biosafety Level 1 (BL1) safety rules.

Safety Notes

CAUTION

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

Required Reagents

Table 1 Required Reagents

Description	Vendor and part number
HaloPlex Target Enrichment System Kit	Select the appropriate kit for your probe design from Table 2
Herculase II Fusion Enzyme with dNTPs (100 mM; 25 mM for each nucleotide), 200 reactions*	Agilent p/n 600677
Nuclease-free Water (not DEPC-treated)	Thermo Fisher Scientific p/n AM9930
AMPure XP Reagent 5 mL 60 mL	Beckman Coulter Genomics p/n A63880 p/n A63881
10 M NaOH, molecular biology grade	Sigma, p/n 72068
2 M acetic acid	Fluka, p/n A8976
10 mM Tris-HCl, pH 8.0 or 10 mM Tris-acetate, pH 8.0	General laboratory supplier
100% Ethanol, molecular biology grade	Sigma-Aldrich p/n E7023
Quant-iT dsDNA BR Assay Kit, for use with the Qubit fluorometer 100 assays, 2-1000 ng 500 assays, 2-1000 ng	Thermo Fisher Scientific p/n Q32850 Thermo Fisher Scientific p/n Q32853

* Also available separately as Herculase II Fusion DNA Polymerase, 40 reactions (Agilent p/n 600675) and 100 mM dNTP Mix (Agilent p/n 200415, sufficient for 1000 enrichment reactions).

Each HaloPlex Target Enrichment System Reagent Kit contains a specific HaloPlex probe. Reagent kit ordering information is provided at [Agilent's SureDesign](#) website and is summarized in [Table 2](#) below.

Table 2 HaloPlex Target Enrichment System Kits for Illumina Sequencing

Probe Design	Part Number	
	96 Reactions	48 Reactions
HaloPlex Custom Panel Tier 1*, ILM	G9901B	G9901C
HaloPlex Custom Panel Tier 1 Plus*, ILM	G9961B	G9961C
HaloPlex Custom Panel Tier 2†, ILM	G9911B	G9911C
HaloPlex Custom Panel Tier 3‡, ILM	G9921B	G9921C

* Tier 1 designs are 1-500 kb with <15,000 probes and require a 3-hour hybridization protocol. Tier 1 Plus designs are also 1-500 kb, but contain 15,000 to 20,000 probes and require a 16-hour hybridization protocol.

† Tier 2 designs are 0.5-2.5 Mb OR 1-500 kb with >20,000 probes.

‡ Tier 3 designs are 2.6 Mb-5 Mb.

NOTE

Kits contain enough reagents for 96 or 48 reactions total, including one or more control reactions using Enrichment Control DNA (ECD) samples. Each run of up to 96 samples should include one ECD control enrichment reaction.

Required Equipment

Table 3 Required Equipment for HaloPlex Target Enrichment

Description	Vendor and part number
Thermal Cycler	Various suppliers *
Thermal cycler-compatible 96-well plates	Consult the thermal cycler manufacturer's recommendations
8-well PCR strip tubes with caps	Agilent p/n 410092 (strip tubes) and Agilent p/n 410096 (strip tube caps)
96-well plate and strip tube-compatible magnetic separator	DynaMag-96 Side magnet, Thermo Fisher Scientific p/n 12331D, or equivalent
1.5 mL tube-compatible magnetic separator	DynaMag-2 magnet, Thermo Fisher Scientific p/n 12321D, or equivalent
Benchtop microcentrifuge	VWR p/n 93000-196, or equivalent
Benchtop plate centrifuge	Labnet International MPS1000 Mini Plate Spinner p/n C1000, or equivalent
Multichannel pipettes (10- μ L and 100- μ L volume)	Pipetman or equivalent
P10, P20, P200 and P1000 pipettes	Pipetman P10, P20, P200, P1000 or equivalent
Adhesive seals for 96-well PCR plates	Agilent p/n 410186, or equivalent
Qubit Fluorometer	Thermo Fisher Scientific p/n Q33226
Qubit assay tubes	Thermo Fisher Scientific p/n Q32856
Ice bucket	General laboratory supplier
Vortex mixer	General laboratory supplier

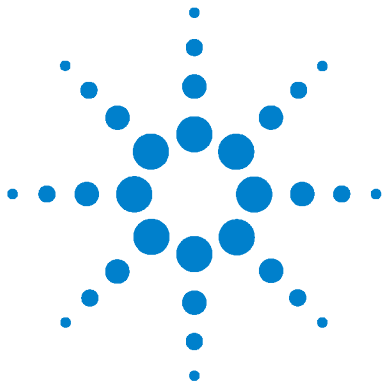
* Thermal cycler must have a maximum reaction volume specification of at least 100 μ L and must be compatible with 0.2 mL tubes.

Optional Validation Reagents and Equipment

Table 4 Reagents and Equipment for Optional Validation Methods

Description	Vendor and part number
Agilent 4200 TapeStation	Agilent p/n G2991AA
96-well sample plates	Agilent p/n 5042-8502
96-well plate foil seals	Agilent p/n 5067-5154
8-well tube strips	Agilent p/n 401428
8-well tube strip caps	Agilent p/n 401425
High Sensitivity D1000 ScreenTape	Agilent p/n 5067-5584
High Sensitivity D1000 Reagents	Agilent p/n 5067-5585
Agilent 2100 Bioanalyzer Instrument	Agilent p/n G2939BA*
High Sensitivity DNA Kit	Agilent p/n 5067-4626
Gel Electrophoresis Platform and Consumables	
XCell SureLock Mini-cell	Thermo Fisher Scientific p/n EI0001
Novex 6% Polyacrylamide, TBE Pre-cast Gels	Thermo Fisher Scientific p/n EC62655BOX
Novex TBE Running Buffer, 5X	Thermo Fisher Scientific p/n LC6675
Novex High-density TBE Sample Buffer, 5X	Thermo Fisher Scientific p/n LC6678
GelRed Nucleic Acid Stain, 3X in water	Biotium p/n 41001
DNA molecular weight markers	General laboratory supplier
UV-transilluminator	General laboratory supplier

* The Agilent 2100 Bioanalyzer Instrument is no longer available for sale. Laboratories equipped with this instrument can purchase and use the associated High Sensitivity DNA Kit for the validation steps detailed in this publication.



2 Sample Preparation

Run Size Considerations	14
Run Time Considerations	14
Step 1. Digest genomic DNA with restriction enzymes	15
Step 2. Hybridize digested DNA to HaloPlex probe for target enrichment and sample indexing	25
Step 3. Capture the target DNA	28
Step 4. Ligate the captured, circularized fragments	31
Step 5. Prepare the PCR Master Mix	32
Step 6. Elute captured DNA with NaOH	33
Step 7. PCR amplify the captured target libraries	34
Step 8. Purify the amplified target libraries	35
Step 9. Validate enrichment and quantify enriched target DNA	37
Step 10. Pool indexed samples and perform multiplexed sequencing	41

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

The target region can vary from 1 kb to 5 Mb. Information on available HaloPlex probe designs can be found at [Agilent's SureDesign](#) website using the *Find Designs* tab.

The HaloPlex Target Enrichment System amplifies thousands of targets in the same reaction, incorporating standard Illumina paired-end sequencing motifs in the process. During hybridization, each sample can be uniquely indexed, allowing for pooling of up to 96 samples per sequencing lane.

See [Figure 1](#) for a summary of the overall HaloPlex target enrichment workflow.



1) Digest genomic DNA.



2) Hybridize the HaloPlex probe library in presence of the Indexing Primer Cassette. Hybridization results in gDNA fragment circularization and incorporation of indexes and Illumina sequencing motifs.



3) Capture target DNA-probe hybrids. Biotinylation of probe DNA allows capture using streptavidin-coated magnetic beads.



4) PCR amplify targeted fragments to produce a sequencing-ready, target-enriched sample.

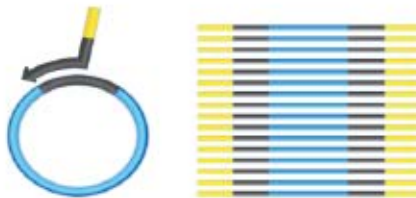


Figure 1 Overall HaloPlex target-enriched sequencing sample preparation workflow.

Run Size Considerations

Kits contain enough reagents for 48 or 96 reactions total, including control reactions using the provided Enrichment Control DNA (ECD). Each run that uses independently-prepared reagent master mixes should include one ECD control enrichment reaction.

The following protocol includes volumes appropriate for 12-sample runs. When planning a run size different from 12 samples, you will need to adjust volumes of components accordingly. Calculate the amount of each solution needed for the number of reactions in your run, plus 2 reactions excess for the restriction digestion steps and 1 reaction excess for the remaining steps. For example, for a 24 reaction run, calculate amounts of each solution by multiplying the single reaction value by 26 for restriction digestion steps and by 25 for hybridization and later steps.

A 96-reaction kit contains enough reagents to prepare master mixes for eight runs of 12 samples each for a total of 96 samples. When processing samples using runs with fewer than 12 samples, some reagents may be depleted before 96 samples are run.

A 48-reaction kit contains enough reagents to prepare master mixes for four runs of 12 samples each, for a total of 48 samples. When processing samples using runs with fewer than 12 samples, some reagents may be depleted before 48 samples are run.

Run Time Considerations

Before you begin, refer to the Certificate of Analysis provided with Box 1 of your kit to determine the hybridization duration appropriate for your design. After reviewing the duration of this and other steps in the protocol, plan the start time for your experiment accordingly.

Designs containing <15,000 probes use a 3-hour hybridization time, and DNA digestion through PCR steps (see [Figure 1](#)) are typically run in the same day. Designs containing $\geq 15,000$ probes use a 16-hour hybridization time, which is typically completed overnight, with the DNA digestion step started in the afternoon.

Step 1. Digest genomic DNA with restriction enzymes

In this step, gDNA samples are digested by 16 different restriction enzymes to create a library of gDNA restriction fragments.

NOTE

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

For HaloPlex target enrichment of FFPE-derived DNA samples, see Agilent publication [G9900-90050](#). This publication provides a PCR-based protocol for assessment of DNA integrity and provides HaloPlex protocol modifications for improved performance from lower-quality DNA samples.

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

Follow the manufacturers instructions for the kits and instruments.

NOTE

Use a fluorometry-based DNA quantitation method, such as Qubit fluorometry or PicoGreen staining, to accurately quantify the DNA starting material.

In the protocol below, 200 ng genomic DNA is split among eight different restriction digests, with an additional 25 ng excess DNA included to allow for pipetting losses. Using <225 ng DNA in the enrichment protocol can result in low yield and can potentiate rare allele dropouts.

- 2 Prepare the DNA samples for the run. For 12-reaction runs, prepare 11 gDNA samples and one Enrichment Control DNA sample.
 - a In separate 0.2-mL PCR tubes, dilute 225 ng of each gDNA sample in 45 μ L nuclease-free water, for a final DNA concentration of 5 ng/ μ L. Store on ice.
 - b In a separate 0.2-mL PCR tube, dispense 45 μ L of the supplied Enrichment Control DNA (ECD). Store on ice.

NOTE

Although specific instructions are provided for the typical 12-sample run size, runs may include up to 96 samples. Include one ECD control sample per run of 2-96 samples. See [page 14](#) for additional run size considerations.

Step 1. Digest genomic DNA with restriction enzymes

3 Prepare the Restriction Enzyme Master Mix strip.

The gDNA is digested in eight different restriction reactions, each containing two restriction enzymes. **The 16 restriction enzymes are provided in two 8-well strip tubes that are distinguished by red and green color markers.** Enzymes are combined from corresponding wells of the red- and green-marked strip tubes, along with restriction buffer and albumin to make eight different RE Master Mixes. [Figure 2](#) illustrates how to prepare the 8-well Restriction Enzyme Master Mix strip for a 12-sample run using the steps detailed on [page 17](#).

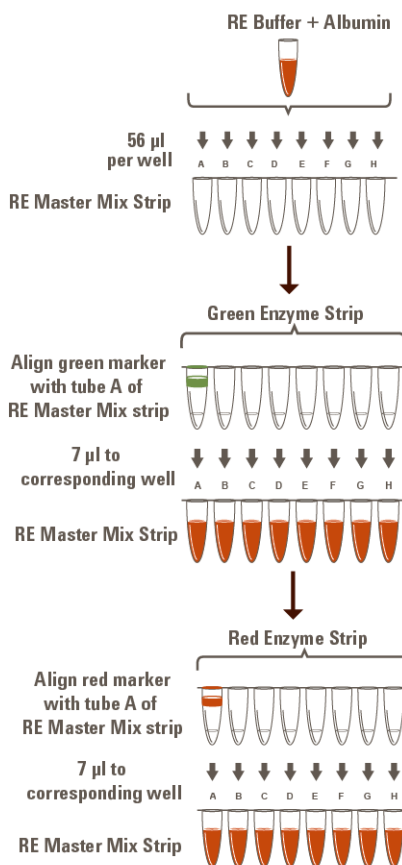


Figure 2 Preparation of the Restriction Enzyme Master Mix Strip for 12-sample run.

Step 1. Digest genomic DNA with restriction enzymes

- a Combine the amounts of RE Buffer and Albumin Solution indicated in the table below in a 1.5-mL tube. Mix by vortexing briefly.

Reagent	Volume for 1 Reaction	Volume for 12 Reactions (includes excess)
RE Buffer	34.0 μ L	476 μ L
re-Albumin Solution*	0.85 μ L	11.9 μ L
Total Volume	34.85 μL	487.9 μL

* re-Albumin Solution replaces the BSA Solution provided with earlier versions of this kit. An equivalent volume of either reagent may be used in this reaction.

- b To begin preparation of the Restriction Enzyme Master Mix Strip, dispense the appropriate volume of the RE Buffer/Albumin mixture to each well of an 8-well strip tube.

Reagent	Volume for 1 Reaction	Volume for 12 Reactions (includes excess)
RE Buffer/Albumin mix	4 μ L	56 μ L

CAUTION

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

- c Using a multichannel pipette, add the appropriate volume of each enzyme from the Green Enzyme Strip, with green marker aligned with tube A, to corresponding tubes A to H of the Restriction Enzyme Master Mix Strip.

Reagent	Volume for 1 Reaction	Volume for 12 Reactions (includes excess)
RE Enzymes from Green Enzyme Strip	0.5 μ L	7 μ L

Step 1. Digest genomic DNA with restriction enzymes

- d** Using a multichannel pipette, add the appropriate volume of each enzyme from the Red Enzyme Strip, with red marker aligned with tube A, to each corresponding tube A to H of the same Restriction Enzyme Master Mix Strip.

Reagent	Volume for 1 Reaction	Volume for 12 Reactions (includes excess)
RE Enzymes from Red Enzyme Strip	0.5 μ L	7 μ L

- e** Mix by gentle vortexing and then spin briefly.
- f** Keep the Restriction Enzyme Master Mix Strip on ice until it is used in [step 4](#).
- 4** Aliquot the Restriction Enzyme Master Mixes to the rows of a 96-well plate to be used as the restriction digest reaction plate.
- a** Align the Restriction Enzyme Master Mix Strip, prepared in [step 3](#), along the vertical side of a 96-well PCR plate as shown below.
- b** Using a multichannel pipette, carefully distribute 5 μ L of each RE master mix row-wise into each well of the plate.

For runs with >12 samples, continue distributing 5 μ L from the same RE Master Mix strip row-wise into each well of the additional plates.

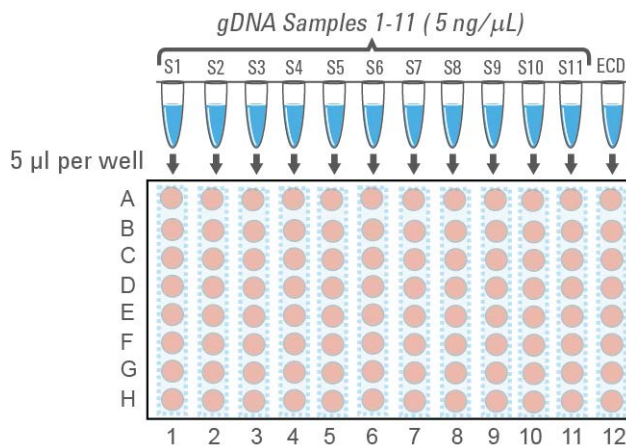
Visually inspect pipette tips for equal volumes before dispensing to the plate(s).



Each row of the 96-well plate now contains 5 μ L per well of the same restriction enzyme combination.

Step 1. Digest genomic DNA with restriction enzymes

- 5** Aliquot DNA samples into the 96-well Restriction Digest Reaction Plate(s).
- a** Align the DNA samples (11 gDNA samples and the ECD sample), prepared in [step 2](#), along the horizontal side of the digestion reaction plate(s) as shown below.



- b** Carefully distribute 5 µL of DNA samples column-wise into each well of the digestion reaction plate.
- If using a multichannel pipette, visually inspect pipette tips for equal volumes before dispensing.
- c** Seal the plate thoroughly with adhesive plastic film.
- 6** Carefully vortex the plate to mix the digestion reactions.
- 7** Briefly spin the plate in a plate centrifuge.

Wells of the prepared 96-well plate now contain complete 10-µL restriction digestion reactions. In this format, each column corresponds to one DNA sample digested in eight different restriction reactions.

Step 1. Digest genomic DNA with restriction enzymes

- 8 Place the Restriction Digest Reaction Plate in a thermal cycler and run the program in [Table 5](#), using a heated lid.

Table 5 Thermal cycler program for restriction digestion

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

NOTE

Do **not** pool the eight restriction digests for a single DNA sample at this time. Restriction enzymes are still active and will catalyze inappropriate cleavage events if DNA samples are pooled before enzyme inactivation. DNA samples are pooled during the hybridization step on [page 25](#), upon which restriction enzymes are inactivated by the reaction conditions.

Step 1. Digest genomic DNA with restriction enzymes

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

Keep the Restriction Digest Reaction Plate on ice during validation.

- a Transfer 4 μ L of each ECD digestion reaction from wells of the digestion reaction plate to fresh 0.2-mL PCR tubes.
- b Incubate the removed 4- μ L samples at 80°C for 5 minutes to inactivate the restriction enzymes.
- c Analyze the prepared samples by electrophoresis using the Agilent 4200 TapeStation (see [page 22](#)), the Agilent 2100 Bioanalyzer (see [page 23](#)), or by gel electrophoresis (see [page 24](#)).

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

NOTE

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

Successful digestion is indicated by the appearance of the three predominant bands. The presence of additional minor bands, with relative abundance similar to the additional bands visible in [Figure 4](#), [Figure 3](#), and [Figure 5](#) does not impact enrichment results.

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

Step 1. Digest genomic DNA with restriction enzymes

Option 1: Validation by Agilent 4200 TapeStation analysis

Use a High Sensitivity D1000 ScreenTape and reagent kit. See the [Agilent High Sensitivity D1000 Assay Quick Guide](#) for assay instructions.

- Prepare an undigested DNA gel control by combining 1 μL of the Enrichment Control DNA solution and 1 μL of nuclease-free water.
- Prepare the TapeStation samples as instructed in the assay Quick Guide. Use 2 μL of each ECD sample diluted with 2 μL of High Sensitivity D1000 sample buffer in separate wells of a tube strip for the analysis.

CAUTION

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

- Load the sample tube strip, the High Sensitivity D1000 ScreenTape, and loading tips into the 4200 TapeStation and start the run. See [Figure 3](#) for sample TapeStation electrophoresis results.

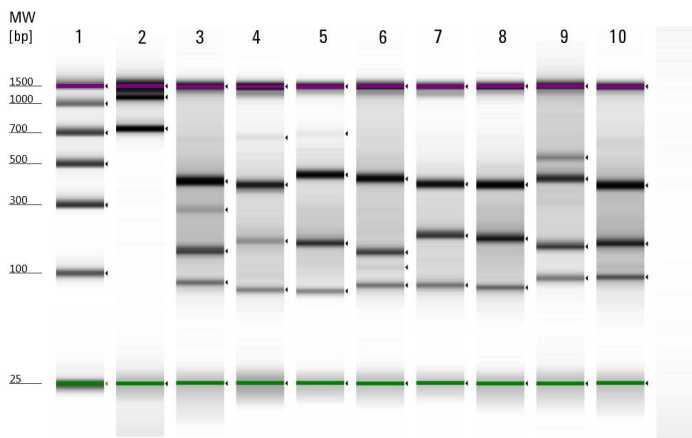


Figure 3 Validation of restriction digestion by TapeStation analysis. Lane 1: High-Sensitivity Ladder, Lane 2: Undigested Enrichment Control DNA, Lanes 3–10: ECD digestion reactions A–H.

Step 1. Digest genomic DNA with restriction enzymes

Option 2: Validation by Agilent 2100 Bioanalyzer analysis

Use a High Sensitivity DNA Kit (p/n 5067-4626) and the 2100 Bioanalyzer system. See the *High Sensitivity DNA Kit Guide* for assay instructions.

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

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

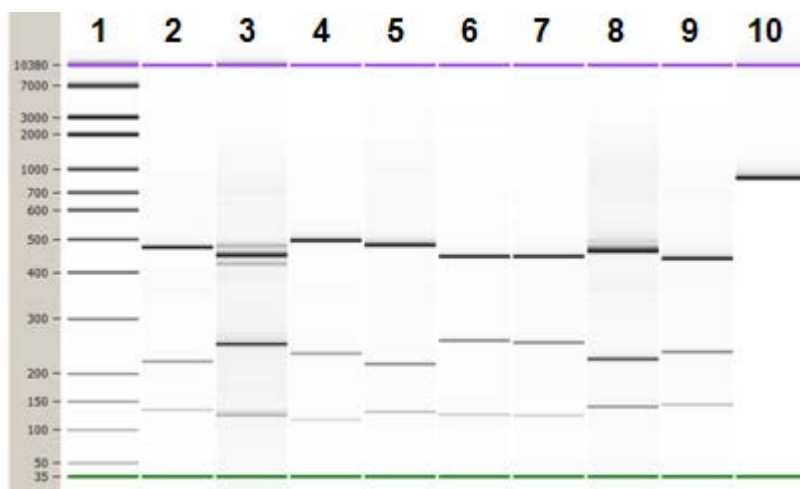


Figure 4 Validation of restriction digestion by 2100 Bioanalyzer system analysis. Lane 1: 50-bp DNA ladder, Lanes 2-9: ECD digestion reactions A–H, Lane 10: Undigested Enrichment Control DNA.

Step 1. Digest genomic DNA with restriction enzymes

Option 3: Validation by gel electrophoresis

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

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

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

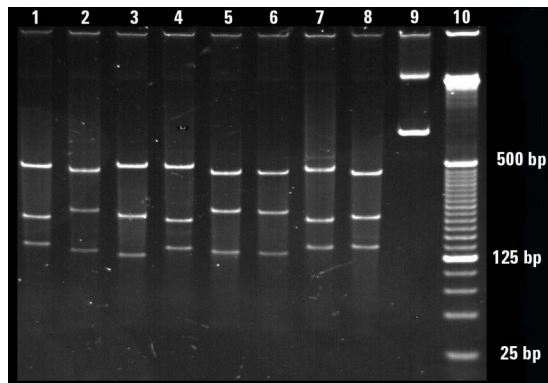


Figure 5 Validation of restriction digestion by gel electrophoresis. Lanes 1–8: ECD digestion reactions A–H, Lane 9: Undigested Enrichment Control DNA, Lane 10: 25-bp DNA ladder.

Stopping Point If you do not continue to the next step, samples may be stored at -20°C for long-term storage. There are no more long-term stopping points until after the PCR amplification step on [page 34](#).

Step 2. Hybridize digested DNA to HaloPlex probe for target enrichment and sample indexing

In this step, the collection of gDNA restriction fragments is hybridized to the HaloPlex probe. The duration of the hybridization reaction is determined by the probe density of your design. Refer to the Certificate of Analysis provided with Box 1 of your kit to determine the hybridization conditions appropriate for your design.

HaloPlex probes are designed to hybridize selectively to fragments originating from target regions of the genome and to direct circularization of the targeted DNA fragments. During the hybridization process, Illumina sequencing motifs including index sequences are incorporated into the targeted fragments.

- 1 Prepare a Hybridization Master Mix by combining the reagents in [Table 6](#). Mix well by gentle vortexing, then spin the tube briefly.

Table 6 Hybridization Master Mix

Reagent	Volume for 1 Reaction	Volume for 12 Reactions (includes excess)
Hybridization Solution	50 μ L	650 μ L
HaloPlex Probe	20 μ L	260 μ L
Total Volume	70 μL	910 μL

- 2 Distribute 70 μ L of the Hybridization Master Mix to each of 12 0.2-mL tubes.
- 3 Add 10 μ L of the appropriate Indexing Primer to each tube containing Hybridization Master Mix.
Be sure to add only one specific Indexing Primer to each hybridization tube, using different indexes for each sample to be multiplexed. Record the identity of the Indexing Primer added to each tube for later sequence analysis.
- 4 Transfer digested DNA samples from the 96-well Restriction Digest Reaction Plate(s) directly into the hybridization reaction tubes prepared in [step 3](#).

Step 2. Hybridize digested DNA to HaloPlex probe for target enrichment and sample indexing

Transfer all eight digestion reactions that correspond to one DNA sample into the appropriate hybridization reaction tube. After addition of each individual digest reaction to the hybridization solution, mix by pipetting before adding the next digest reaction to ensure complete inactivation of the enzymes.

CAUTION

Do **not** pool the digestion samples before adding to the hybridization reaction mixture as restriction enzymes are still active and may catalyze inappropriate cleavage events.

For the ECD sample, add 32 μL of nuclease-free water, in addition to the digested DNA samples, to compensate for the volume removed for digest validation.

After pooling, each hybridization reaction contains the following components:

- 70 μL Hybridization Master Mix
- 10 μL Indexing Primer
- approximately 80 μL pooled digested DNA samples

NOTE

Due to partial evaporation of samples, you may recover less than 10 μL of each restriction digest. Minor reductions to the digested DNA pool volume will not impact hybridization performance; you do not need to compensate for any sample evaporation volume losses in the final pool.

5 Vortex the mixtures briefly and then spin tubes briefly.

Step 2. Hybridize digested DNA to HaloPlex probe for target enrichment and sample indexing

- 6 Place the hybridization reaction tubes in a thermal cycler. Run the appropriate program in Table 7, using the hybridization duration listed on the Certificate of Analysis.

Use a heated lid. Do **not** include a low-temperature hold step in the thermal cycler program. Incubation at 54°C for more than the indicated time is not recommended.

Table 7 Thermal cycler program* for probe hybridization

Step	Temperature	Time (Duration of Step)	
		Designs with <15,000 probes (see Certificate of Analysis)	Designs with ≥15,000 probes (see Certificate of Analysis)
Step 1	95°C	10 minutes	10 minutes
Step 2	54°C	3 hours	16 hours

* Thermal cyclers that use calculated temperature methods cannot be set to 160 µL reaction volumes. In that case, enter the maximum possible volume.

CAUTION

Make sure that the thermal cycler has a maximum reaction volume specification of at least 100 µL. The performance of the selected thermal cycler in this application should be verified before use.

Step 3. Capture the target DNA

In this step, the circularized target DNA-probe hybrids, containing biotin, are captured on streptavidin beads.

- 1 Remove reagents to be used in upcoming protocol steps from cold storage and allow the solutions to reach room temperature:
 - From -20°C storage, remove the Capture Solution, Wash Solution, Ligation Solution and SSC Buffer.
 - From $+4^{\circ}\text{C}$ storage, remove the HaloPlex Magnetic Beads.
- 2 Obtain or prepare 0.5 μL per sample, plus excess, of 2 M acetic acid, for use on [page 32](#).

NOTE

Use a high-quality 2 M acetic acid solution. See [Table 1](#) on page 8 for 2 M acetic acid supplier information, or prepare 2 M acetic acid from high-quality glacial acetic acid.

- 3 Prepare 25 μL per sample, plus excess, of fresh 50 mM NaOH for use in the DNA elution step on [page 33](#).

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

CAUTION

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

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

- 4 Vigorously resuspend the provided HaloPlex Magnetic Beads on a vortex mixer. The magnetic beads settle during storage.

5 Prepare 40 μL (1 Volume) of HaloPlex Magnetic Beads per hybridization sample, plus excess, for the capture reaction:

a Transfer the appropriate volume of bead suspension to a 1.5-mL tube.

Reagent	Volume for 1 Reaction	Volume for 12 Reactions (includes excess)
HaloPlex Magnetic Bead suspension	40 μL	520 μL

b Put the tube into a 1.5 mL tube-compatible magnetic rack for 5 minutes.

c After verifying that the solution has cleared, carefully remove and discard the supernatant using a pipette.

d Add an equivalent volume of Capture Solution to the beads and resuspend by pipetting up and down.

Reagent	Volume for 1 Reaction	Volume for 12 Reactions (includes excess)
Capture Solution	40 μL	520 μL

6 Remove the hybridization reactions from the thermal cycler and immediately add 40 μL of the prepared bead suspension to each 160- μL hybridization reaction.

NOTE

When adding beads to the hybridization reactions, visually inspect the bead preparation to ensure a homogeneous suspension with no aggregated bead mass at the bottom of the tube. If aggregation is present, thoroughly resuspend the beads by vortexing and pipetting up and down before use.

7 After adding the magnetic beads, mix the capture reactions thoroughly by pipetting up and down 15 times using a 100- μL pipette set to 80 μL .

8 Incubate the capture reactions at room temperature for 15 minutes.

9 Briefly spin the tubes in a desktop centrifuge and then transfer the tubes to a strip tube-compatible magnetic plate.

NOTE

Use the magnetic plate for the remainder of magnetic bead collection steps for samples in PCR tubes or strip tubes.

- 10** Wait for the solution to clear (about 30 seconds), then remove and discard the supernatant using a pipette set to 200 μL .
- 11** Wash the bead-bound samples:
 - a** Remove the capture reaction tubes from the magnetic plate and add 100 μL of Wash Solution to each tube.
 - b** Resuspend the beads thoroughly by pipetting up and down 10 times using a 100- μL multichannel pipette set to 80 μL .
 - c** Incubate the tubes in a thermal cycler at 46°C for 10 minutes, using a heated lid.

Do **not** include a low-temperature hold step in the thermal cycler program following the 10-minute incubation.
 - d** Briefly spin the tubes in a desktop centrifuge at room temperature and then transfer the tubes to the magnetic plate.
 - e** Wait for the solution to clear (about 30 seconds), then carefully remove and discard the supernatant using a pipette set to 120 μL . If necessary, carefully remove any residual liquid with a 20- μL volume pipette.

Step 4. Ligate the captured, circularized fragments

In this step, DNA ligase is added to the capture reaction to close nicks in the circularized probe-target DNA hybrids.

- 1 Prepare a DNA ligation master mix by combining the reagents in the following table.

Mix the components thoroughly by gentle vortexing then spin the tube briefly.

Table 8 Preparation of DNA ligation master mix

Reagent	Volume for 1 Reaction	Volume for 12 Reactions (includes excess)
Ligation Solution	47.5 μ L	617.5 μ L
DNA Ligase	2.5 μ L	32.5 μ L
Total Volume	50 μL	650 μL

- 2 Add 50 μ L of the DNA ligation master mix to the beads in each DNA capture reaction tube.
- 3 Resuspend the beads thoroughly by pipetting up and down 15 times using a 100- μ L multichannel pipette set to 40 μ L.
- 4 Incubate the tubes in a thermal cycler at 55°C for 10 minutes, using a heated lid.

The thermal cycler may be programmed to include a 4°C hold step following the 10-minute incubation.

During the 10-minute incubation, prepare the PCR master mix as specified in the following step.

Step 5. Prepare the PCR Master Mix

In this step, you prepare a PCR master mix for the captured target DNA amplification step on [page 34](#).

CAUTION

It is critical to include Acetic acid at 2 M concentration in this step to ensure neutralization of the NaOH used for elution on [page 33](#).

- 1 Prepare the PCR master mix by combining the reagents in the following table.

Table 9 Preparation of PCR master mix

Reagent	Volume for 1 reaction	Volume for 12 reactions (includes excess)
Nuclease-free water	16.1 μL	209.3 μL
5X Herculase II Reaction Buffer	10 μL	130 μL
dNTPs (100 mM, 25 mM for each dNTP)	0.4 μL	5.2 μL
Primer 1 (25 μM)	1 μL	13 μL
Primer 2 (25 μM)	1 μL	13 μL
2 M Acetic acid	0.5 μL	6.5 μL
Herculase II Fusion DNA Polymerase	1 μL	13 μL
Total	30 μL	390 μL

- 2 Mix the master mix components by gentle vortexing, then distribute 30- μL aliquots to fresh 0.2-mL reaction tubes.
- 3 Store the tubes on ice until they are used in “[Step 7. PCR amplify the captured target libraries](#)” on [page 34](#).

Step 6. Elute captured DNA with NaOH

When the 10-minute ligation reaction period is complete, proceed with the following steps to elute the captured DNA libraries.

CAUTION

Using a high-quality NaOH solution for this step is critical for optimal DNA elution and recovery.

Be sure to use freshly-prepared 50 mM NaOH, prepared from 10 M NaOH according to the instructions on [page 28](#).

- 1 Briefly spin the ligation reaction tubes in a desktop centrifuge and then transfer the tubes to the magnetic plate.
- 2 Wait for the solution to clear (about 30 seconds), then carefully remove and discard the supernatant using a pipette set to 50 μ L.
- 3 Remove the tubes from the magnetic plate and add 100 μ L of the SSC Buffer provided with the kit to each tube.
- 4 Resuspend the beads thoroughly by pipetting up and down 10 times using a 100- μ L multichannel pipette set to 80 μ L.
- 5 Briefly spin the tubes and then return the tubes to the magnetic plate.
- 6 Wait for the solution to clear (about 30 seconds), then carefully remove and discard the SSC Buffer using a multichannel pipette set to 120 μ L. If necessary, carefully remove any residual liquid with a 20- μ L volume pipette.
- 7 Add 25 μ L of 50 mM NaOH, which was freshly-prepared on [page 28](#), to each tube.
- 8 Resuspend the beads thoroughly by pipetting up and down 10 times using a 100- μ L multichannel pipette set to 15 μ L.
- 9 Incubate samples for 1 minute at room temperature to allow elution of the captured DNA.
- 10 Briefly spin the tubes and then transfer the tubes to the magnetic plate. Proceed immediately to PCR amplification in the following section.

Step 7. PCR amplify the captured target libraries

- 1 Prepare amplification reactions by transferring 20 μ L of cleared supernatant from each tube on the magnetic plate to a PCR Master Mix tube held on ice (from [page 32](#)).
- 2 Mix by gentle vortexing and then spin briefly to collect the liquid.
- 3 Place the amplification reaction tubes in a thermal cycler and run the program in [Table 10](#), using a heated lid.

The optimal amplification cycle number varies for each HaloPlex Probe design. Consult the Certificate of Analysis (provided with Box 1 of your kit) for the PCR cycling recommendation for your probe.

Table 10 Post-capture DNA amplification PCR program

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

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

Step 8. Purify the amplified target libraries

In this step, the amplified target DNA is purified using AMPure XP beads.

- 1 Let the AMPure XP beads come to room temperature for at least 30 minutes.
- 2 Prepare 400 μL of 70% ethanol per sample, plus excess, for use in [step 10](#).
- 3 Transfer 40 μL of each PCR reaction sample to a fresh 0.2-mL tube. Store the remaining volume of each sample at -20°C for troubleshooting.
- 4 Mix the AMPure XP bead suspension well, until the suspension appears homogeneous and consistent in color.
- 5 For each sample to be purified, prepare a bead mix by combining 40 μL of nuclease-free water and 100 μL of the homogenous AMPure XP bead suspension. Mix well, until the bead mix suspension appears homogeneous.
- 6 Add 140 μL of the homogeneous bead suspension prepared in [step 5](#) to each 40- μL amplified library sample. Vortex thoroughly.
Using this bead-to-sample volume ratio is imperative to ensure optimal purification results.
- 7 Incubate samples for 5 minutes at room temperature with continuous shaking.
Make sure the samples are properly mixing in the wells during the 5-minute incubation.
- 8 Spin briefly to collect the liquid, then place the tubes in the magnetic plate. Wait for the solution to clear (approximately 5 minutes).
- 9 Keep the tubes in the magnetic plate. Carefully remove and discard the cleared solution from each tube using a 200- μL pipette set to 180 μL . Do not touch the beads while removing the solution.
- 10 Continue to keep the tubes in the magnetic plate while you add 200 μL of 70% ethanol into the tubes.
Use fresh 70% ethanol for optimal results.
- 11 Wait for 30 seconds to allow any disturbed beads to settle, then remove the ethanol using a 200- μL pipette set to 200 μL .
- 12 Repeat [step 10](#) and [step 11](#) once for a total of two washes.
- 13 Remove any residual ethanol with a 20- μL volume pipette.

14 Air-dry the tubes with open lids at room temperature until the residual ethanol completely evaporates.

Make sure all ethanol has evaporated before continuing.

15 Remove tubes from the magnetic plate and add 40 μL of 10 mM Tris-acetate or Tris-HCl buffer (pH 8.0) to each sample.

NOTE

Use room-temperature Tris-acetate or Tris-HCl buffer for elution at this step.

16 Mix thoroughly by pipetting up and down 15 times using a 100- μL pipette set to 30 μL .

17 Incubate for 2 minutes at room temperature to allow elution of DNA.

18 Put the tube in the magnetic plate and leave for 2 minutes or until the solution is clear.

19 Remove the cleared supernatant (approximately 40 μL) to a fresh tube. You can discard the beads at this time.

Stopping Point

If you do not continue to the next step, samples may be stored at -20°C for long-term storage (up to one year). Avoid subjecting the stored DNA samples to multiple freeze-thaw cycles.

Step 9. Validate enrichment and quantify enriched target DNA

Prior to sample pooling and sequencing sample preparation, validate enrichment and quantify the enriched target DNA in each library sample using the Agilent 4200 TapeStation (see [page 38](#)) or 2100 Bioanalyzer (see [page 38](#)).

Enriched library samples may also be qualitatively analyzed using gel electrophoresis. Sample gel electrophoresis results are provided in the [Reference](#) section on [page 47](#).

Expected Results

Each amplicon in the prepared library contains one target insert surrounded by sequence motifs required for multiplexed sequencing using the Illumina platform. Amplicons include 50 to 500 bp of target DNA insert and 125 bp of sequencing motifs, as shown in [Figure 6](#).



Figure 6 Content of target-enriched amplicons. Each amplicon contains one target insert (blue) surrounded by the Illumina paired-end sequencing elements (black), the sample index (red) and the library bridge PCR primers (yellow).

The amplicons should range from 175 to 625 bp in length, with the majority of products sized 225 to 525 bp. Amplicons in the 175 to 625 bp size range should be included for quantitation of the enriched target DNA in each sample. Any spurious DNA products outside of this size range in any sample should be excluded from the target DNA quantitation results.

Step 9. Validate enrichment and quantify enriched target DNA

Option 1: Analysis using the Agilent 4200 TapeStation

Use a High Sensitivity D1000 ScreenTape and reagent kit. See the [Agilent High Sensitivity D1000 Assay Quick Guide](#) for assay instructions.

- 1 Prepare the TapeStation samples as instructed in the reagent Quick Guide. Use 2 μL of each enriched library sample diluted with 2 μL of High Sensitivity D1000 sample buffer in separate wells of a tube strip for the analysis.

CAUTION

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

- 2 Load the sample tube strip, the High Sensitivity D1000 ScreenTape, and loading tips into the 4200 TapeStation and start the run.
- 3 Analyze the electropherogram for each sample using the analysis guidelines on [page 40](#).

See [Figure 7](#) for a sample electropherogram.

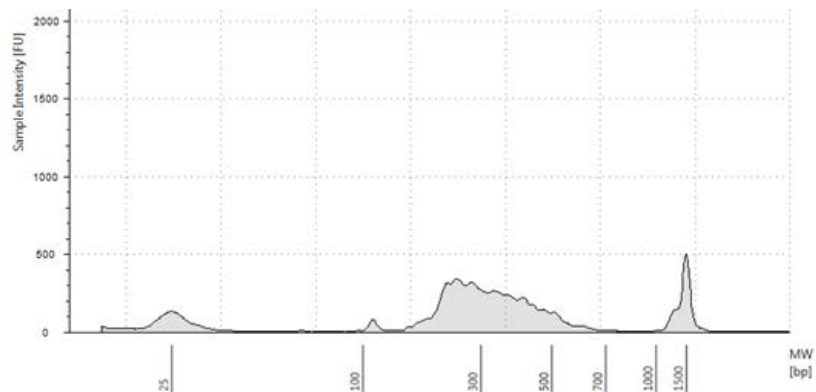


Figure 7 Validation of enrichment by TapeStation analysis.

Step 9. Validate enrichment and quantify enriched target DNA

Option 2: Analysis using the Agilent 2100 Bioanalyzer

Use a Bioanalyzer High Sensitivity DNA Assay kit and the Agilent 2100 Bioanalyzer. See the *High Sensitivity DNA Kit Guide* for assay instructions.

- 1 Prepare the chip, samples and ladder as instructed in the reagent kit guide, using 1 μL of enriched library sample for the analysis.
- 2 Load the prepared chip into the instrument and start the run within five minutes after preparation.
- 3 Analyze the electropherogram for each sample using the analysis guidelines on [page 40](#).

See [Figure 8](#) for a sample electropherogram.

NOTE

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

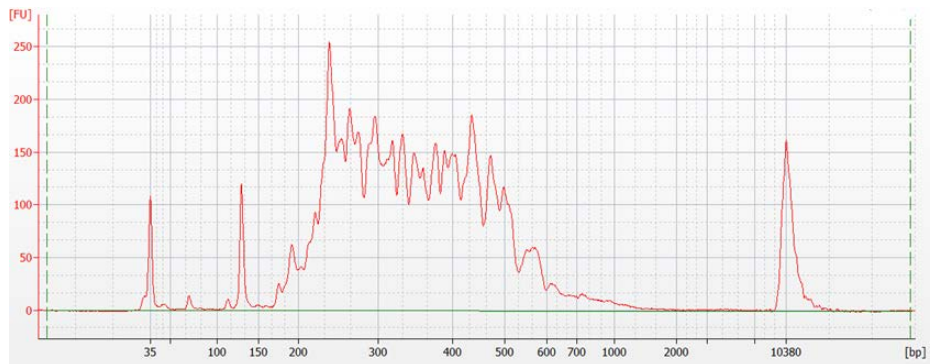


Figure 8 Validation of enrichment by Agilent 2100 Bioanalyzer analysis.

Step 9. Validate enrichment and quantify enriched target DNA

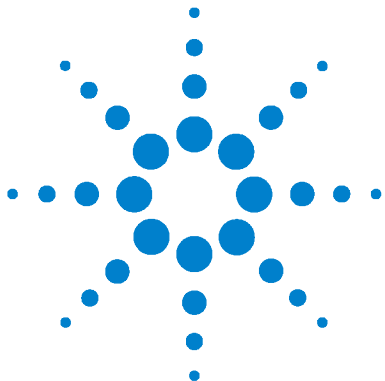
Analysis of Electropherogram Results

- Check that the electropherogram shows a peak fragment size between approximately 225 to 525 bp.
- Determine the concentration of enriched target DNA in the sample by integration under the peak between 175 and 625 bp. Peaks at <150 bp may be observed, but should be excluded from quantitation.
- Some designs may generate a peak at about 125 bp. This peak is associated with an adaptor-dimer product which will cluster and generate sequence that does not map to the genome. If the molar fraction of the 125 bp peak is greater than 10%, do another round of AMPure purification after pooling samples. First, pool equimolar amounts of libraries to be multiplexed, using concentrations determined for the 175–625 peak of each sample. Using 40 μ L of the pooled libraries, purify the DNA using AMPure XP beads according to the protocol on [page 35](#).

Step 10. Pool indexed samples and perform multiplexed sequencing

Use the following guidelines to design your sample pooling and sequencing strategy:

- Use the TapeStation- or Bioanalyzer-measured concentration of 175-625 bp products in each sample to pool equimolar amounts of differentially indexed samples in order to optimize the use of sequencing capacity.
- The final HaloPlex enrichment pool is ready for direct sequencing using standard Illumina paired-end primers and chemistry on an Illumina platform. If you need assistance with the NGS segment of the workflow, please contact Agilent's NGS Support team or your local representative.
- Use 100 + 100 bp or 150 + 150 bp paired-end sequencing, depending on the selection made during probe design. Since the read length affects maximum achievable coverage, check the design report to verify read length selected in probe design.
- Sequencing runs must be set up to perform an 8-nt Index 1 (i7) read.
- For complete HaloPlex index sequence information see [page 46](#).
- Before aligning reads to the reference genome, trim the Illumina adaptor sequences from the reads.



3 Reference

Kit Contents 43
Nucleotide Sequences of HaloPlex Indexes 46
Qualitative analysis of enrichment by gel electrophoresis 47

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



Kit Contents

HaloPlex Target Enrichment System Kits for custom designs include the component kits listed in [Table 11](#).

Table 11 HaloPlex Target Enrichment System Kit Contents-Custom Designs

Design Type	Kit Part Number	HaloPlex Target Enrichment System-ILM, Box 1*	HaloPlex Magnetic Beads Box 2
		Store at -20°C	Store at +4°C
HaloPlex 1-500 kb with <15,000 Probes, ILM	G9901C (48 Reactions)	5190-8044 OR 5190-8045	5190-5976
	G9901B (96 Reactions)	5190-8050 OR 5190-8051	5190-5386
HaloPlex 1-500 kb with 15,000-20,000 Probes, ILM	G9961C (48 Reactions)	5191-4066 OR 5191-4067	5190-5976
	G9961B (96 Reactions)	5191-4068 OR 5191-4069	5190-5386
HaloPlex 0.5-2.5 Mb OR <0.5 Mb with >20,000 probes, ILM	G9911C (48 Reactions)	5190-8046 OR 5190-8047	5190-5976
	G9911B (96 Reactions)	5190-8052 OR 5190-8053	5190-5386
HaloPlex 2.6 Mb-5 Mb, ILM	G9921C (48 Reactions)	5190-8048 OR 5190-8049	5190-5976
	G9921B (96 Reactions)	5190-8054 OR 5190-8055	5190-5386

* See [Table 12](#) for list of included reagents. Part number 5190-8044, 5190-8050, 5191-4066, 5191-4068, 5190-8046, 5190-8052, 5190-8048 or 5190-8054 is provided for the first order of a specific HaloPlex Custom Probe design. Re-order kits, containing previously-purchased Custom Probe designs, include Box 1 part number 5190-8045, 5190-8051, 5191-4067, 5191-4069, 5190-8047, 5190-8053, 5190-8049 or 5190-8055.

The contents of the HaloPlex Target Enrichment System Box 1 included with each kit are detailed in [Table 12](#).

Table 12 HaloPlex Target Enrichment System Box 1 Contents

Included Reagents	48 Reaction Kit	96 Reaction Kit
Hybridization Solution	bottle	bottle
Ligation Solution	bottle	bottle
Wash Solution	bottle	bottle
Capture Solution	bottle	bottle
SSC Buffer	bottle	bottle
RE Buffer	bottle	bottle
re-Albumin Solution*	tube with clear cap	tube with clear cap
DNA Ligase	tube with red cap	tube with red cap
Enrichment Control DNA	tube with orange cap	tube with orange cap
Primer 1	tube with yellow cap	tube with yellow cap
Primer 2	tube with blue cap	tube with blue cap
HaloPlex Indexing Primers	96-well plate with Indexing Primer A01 to H06 (blue plate) [†]	96-well plate with Indexing Primer A01 to H12 (blue plate) [‡]
Enzyme Strip 1	8-well strip tube with green label	8-well strip tube with green label
Enzyme Strip 2	8-well strip tube with red label	8-well strip tube with red label
HaloPlex Probe	tube with pink cap	tube with pink cap

* re-Albumin Solution replaces the BSA Solution provided with earlier versions of this kit. Either reagent solution can be used in the protocols provided in this publication (see [page 16](#) to [page 17](#)).

† See [Table 13](#) for a plate map.

‡ See [Table 14](#) for a plate map.

Table 13 Plate map for HaloPlex Indexing Primers A01 through H06 provided with 48-reaction kits; wells in columns 7 through 12 are empty

	1	2	3	4	5	6	7	8	9	10	11	12
A	A01	A02	A03	A04	A05	A06	–	–	–	–	–	–
B	B01	B02	B03	B04	B05	B06	–	–	–	–	–	–
C	C01	C02	C03	C04	C05	C06	–	–	–	–	–	–
D	D01	D02	D03	D04	D05	D06	–	–	–	–	–	–
E	E01	E02	E03	E04	E05	E06	–	–	–	–	–	–
F	F01	F02	F03	F04	F05	F06	–	–	–	–	–	–
G	G01	G02	G03	G04	G05	G06	–	–	–	–	–	–
H	H01	H02	H03	H04	H05	H06	–	–	–	–	–	–

Table 14 Plate map for HaloPlex Indexing Primers A01 through H12 provided with 96-reaction kits

	1	2	3	4	5	6	7	8	9	10	11	12
A	A01	A02	A03	A04	A05	A06	A07	A08	A09	A10	A11	A12
B	B01	B02	B03	B04	B05	B06	B07	B08	B09	B10	B11	B12
C	C01	C02	C03	C04	C05	C06	C07	C08	C09	C10	C11	C12
D	D01	D02	D03	D04	D05	D06	D07	D08	D09	D10	D11	D12
E	E01	E02	E03	E04	E05	E06	E07	E08	E09	E10	E11	E12
F	F01	F02	F03	F04	F05	F06	F07	F08	F09	F10	F11	F12
G	G01	G02	G03	G04	G05	G06	G07	G08	G09	G10	G11	G12
H	H01	H02	H03	H04	H05	H06	H07	H08	H09	H10	H11	H12

Nucleotide Sequences of HaloPlex Indexes

The nucleotide sequence of the 8-nucleotide index portion of each HaloPlex Indexing Primer is provided in the table below.

Table 15 HaloPlex Indexes, for indexing primers provided in blue 96-well plate

Index	Sequence	Index	Sequence	Index	Sequence	Index	Sequence
A01	ATGCCTAA	A04	AACTCACC	A07	ACGTATCA	A10	AATGTTGC
B01	GAATCTGA	B04	GCTAACGA	B07	GTCTGTCA	B10	TGAAGAGA
C01	AACGTGAT	C04	CAGATCTG	C07	CTAAGGTC	C10	AGATCGCA
D01	CACTTCGA	D04	ATCCTGTA	D07	CGACACAC	D10	AAGAGATC
E01	GCCAAGAC	E04	CTGTAGCC	E07	CCGTGAGA	E10	CAACCACA
F01	GACTAGTA	F04	GCTCGGTA	F07	GTGTTCTA	F10	TGGAACAA
G01	ATTGGCTC	G04	ACACGACC	G07	CAATGGAA	G10	CCTCTATC
H01	GATGAATC	H04	AGTCACTA	H07	AGCACCTC	H10	ACAGATTC
A02	AGCAGGAA	A05	AACGCTTA	A08	CAGCGTTA	A11	CCAGTTCA
B02	GAGCTGAA	B05	GGAGAACA	B08	TAGGATGA	B11	TGGCTTCA
C02	AAACATCG	C05	CATCAAGT	C08	AGTGGTCA	C11	CGACTGGA
D02	GAGTTAGC	D05	AAGGTACA	D08	ACAGCAGA	D11	CAAGACTA
E02	CGAACTTA	E05	CGCTGATC	E08	CATACCAA	E11	CCTCCTGA
F02	GATAGACA	F05	GGTGCGAA	F08	TATCAGCA	F11	TGGTGGTA
G02	AAGGACAC	G05	CCTAATCC	G08	ATAGCGAC	G11	AACAACCA
H02	GACAGTGC	H05	CTGAGCCA	H08	ACGCTCGA	H11	AATCCGTC
A03	ATCATTCC	A06	AGCCATGC	A09	CTCAATGA	A12	CAAGGAGC
B03	GCCACATA	B06	GTACGCAA	B09	TCCGTCTA	B12	TTCACGCA
C03	ACCACTGT	C06	AGTACAAG	C09	AGGCTAAC	C12	CACCTTAC
D03	CTGGCATA	D06	ACATTGGC	D09	CCATCCTC	D12	AAGACGGA
E03	ACCTCCAA	E06	ATTGAGGA	E09	AGATGTAC	E12	ACACAGAA
F03	GCGAGTAA	F06	GTCGTAGA	F09	TCTTCACA	F12	GAACAGGC
G03	ACTATGCA	G06	AGAGTCAA	G09	CCGAAGTA	G12	AACCGAGA
H03	CGGATTGC	H06	CCGACAAC	H09	CGCATACA	H12	ACAAGCTA

Qualitative analysis of enrichment by gel electrophoresis

Enrichment products may be qualitatively analyzed by gel electrophoresis. Analyze 5 μ L of each enriched library sample (enriched ECD sample or experimental enriched libraries) by electrophoresis on a Novex 6% polyacrylamide TBE pre-cast gel. See [page 21](#) for additional gel analysis protocol recommendations.

Successful enrichment is indicated by the presence of a smear of amplicons from approximately 225 to 525 bp in each enrichment library lane. For some probe designs, low molecular weight (<150 bp) bands may also be visible, but should not be included in enriched sample quantitation. See [Figure 9](#) for a sample gel analysis image.

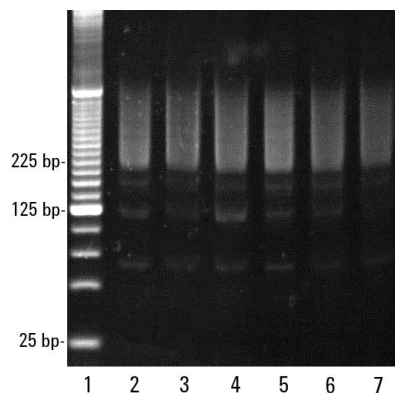


Figure 9 Validation of HaloPlex enrichment process by gel electrophoresis. Lane 1: 25-bp DNA ladder, Lanes 2-7: enriched library samples.

In This Book

This guide contains information to run the HaloPlex Target Enrichment System protocol for the Illumina sequencing platform.

© Agilent Technologies, Inc. 2015, 2018,
2025

Version H0, February 2025



G9900-90001