

Avida Methyl Reagent Kits

For targeted methylation sequencing on the Illumina platform

Library preparation, target capture, conversion, and indexing

Protocol

Version C1, November 2025

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

Notices

© Agilent Technologies, Inc. 2024-2025

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

G9419-90000

Edition

Version C1, November 2025

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 provides an optimized protocol for preparation of target-enriched Illumina paired-end sequencing libraries using Avida Methyl reagent kits.

1 Before You Begin

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

2 gDNA Fragmentation

This chapter describes the steps to fragment gDNA samples using mechanical shearing. gDNA samples must be fragmented prior to starting the Avida Methyl workflow.

3 PCR-Free Library Preparation

This chapter describes the steps for the PCR-free library preparation section of the Avida Methyl workflow. NGS libraries are prepared using adaptors that include Unique Molecular Identifiers (UMIs).

4 Target Capture

This chapter describes the steps for the target capture section of the Avida Methyl workflow. The adaptor-ligated libraries are hybridized with a target-specific panel. After hybridization, the targeted molecules are captured on Capture Beads coated with Streptavidin.

5 Soft Conversion and Repair

This chapter describes the steps for the soft conversion and repair section of the Avida Methyl workflow. The adaptor-ligated libraries are treated with soft conversion reagents to deaminate any unmethylated cytosines, thereby converting them to uracils. The converted DNA is then purified, washed, and repaired. During the PCR indexing step (Chapter 6, "Indexing PCR, Library Purification, and Quality Assessment") the uracils are replaced with thymines in the PCR products.

6 Indexing PCR, Library Purification, and Quality Assessment

This chapter describes the steps for the indexing PCR, library purification, and quality assessment section of the Avida Methyl workflow. The captured libraries are amplified and indexed, and the indexed libraries are then purified with AMPure XP Beads. Finally, the purified, indexed libraries are quantified and analyzed for quality.

7 Sequencing and NGS Analysis

This chapter contains guidance on library sequencing and analysis. Refer to your specific Illumina sequencer's user guide for specific instructions on how to perform sequencing.

8 Reference

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

What's New in Version C1

- Updated information on Avida panel sizes. See [Table 3](#) on page 13 and the accompanying *Note* below the table.
- Simplified preparation of the Library Binding Beads that eliminates 5-minute incubation of the stock tube/bottle prior to use. See [step 2](#) on [page 25](#).
- Instruction to reheat the Hyb Buffer prior to use if a precipitate forms during its storage at room temperature. See [step 2](#) on [page 28](#).
- Increased excess volumes provided in [Table 24](#) on page 36 for preparation of 0.1X Soft Conversion Elution Buffer.
- Additional sequencing run guidelines, including Illumina Kit Configuration Selection Guidelines, added to "[Step 3. Sequence the libraries](#)" on page 49.
- Updated kit configuration information for Soft Conversion Reagent B, which may be shipped as a separate part number or included in the Avida Methyl and Duo Reagent Box 3. See [Table 2](#) on page 12 and [Table 38](#) on page 55.

What's New in Version B2

- Additional information on Avida panel size values and calculations. See [Table 3](#) on page 13, including the accompanying *Note*, for details.

What's New in Version B1

- Corrected kit part numbers in [Table 38](#) on page 55.

What's New in Version B0

- Throughout Protocol, using the term *soft conversion* to refer to the Avida soft conversion process for deamination of unmethylated cytosines.
- Expanded information on compatible reagent kits for the Agilent Fragment Analyzer system, including addition of electropherogram images in [Figure 4](#) and [Figure 5](#).
- In [Quick Reference Protocol](#), clarified storage conditions for the Avida panels and the final sequencing libraries.
- Added *Note* ([page 25](#)) on preparing aliquots of Hyb Mix 1.
- Updated thawing conditions for Adaptor for ILM ([page 20](#) and [page 22](#)).
- Expanded recommended Qubit Assay Kits to include Qubit dsDNA HS and BR Assay Kits.
- Updated the instructions for "[Step 2. Capture Bead Preparation](#)" on page 29, including adding details based on number of reactions.
- Updated storage conditions for the Stopping Point on [page 42](#).
- In the [Quick Reference Protocol](#), added link to the new graphics-based Avida Methyl Quick Start Protocol.

Content

1 Before You Begin

Introduction	8
Avida Target Enrichment Technology	8
Overview of the Workflow	9
Procedural and Safety Notes	10
Important practices for preventing contamination	10
Safety notes	10
Procedural notes	11
Materials Provided with Avida Methyl Kits and Panels	12
Online resource for Avida index sequences	13
Additional Materials Required	14
Input DNA samples	14
Additional reagents and equipment required	14
Equipment required for gDNA fragmentation	16

2 gDNA Fragmentation

Step 1. Prepare for shearing	18
Step 2. Shear the gDNA samples	18

3 PCR-Free Library Preparation

Step 1. End prep (end repair and dA-tailing)	21
Step 2. Ligation	23
Step 3. Library bead binding	25

4 Target Capture

Step 1. Hybridization	28
Step 2. Capture Bead Preparation	29
Step 3. Bead Capture	31
Step 4. Washing (Bead-Bound DNA Hybrids)	32

5 Soft Conversion and Repair

Step 1. Soft Conversion	34
Step 2. Bead Purification	36
Step 3. Washing (Bead-Bound Converted Libraries)	37

Step 4. DNA Repair	38
6 Indexing PCR, Library Purification, and Quality Assessment	
Step 1. Indexing PCR	40
Step 2. Library Purification	42
Step 3. Library Quality Assessment and Quantification	43
7 Sequencing and NGS Analysis	
Step 1. Pool samples for multiplexed sequencing	47
Step 2. Prepare the sequencing samples	48
Step 3. Sequence the libraries	49
Step 4. Process and analyze the reads	51
8 Reference	
Reagent Kit Contents	54
Avida Sub-Kit Details	54
Avida Index Primer Pair Information	56
Index Primer Pair Strip Tube and Plate Maps	56
Avida Custom AddOn Panels	58
Troubleshooting Guide	59
Quick Reference Protocol	60

1

Before You Begin

Introduction [8](#)

Overview of the Workflow [9](#)

Procedural and Safety Notes [10](#)

Materials Provided with Avida Methyl Kits and Panels [12](#)

Additional Materials Required [14](#)

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

NOTE

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

Introduction

DNA methylation changes underlie the mechanisms of many physiological and pathological conditions, including the development of cancer. There is a growing demand in research and clinical settings for a simple, ultra-sensitive targeted sequencing tool for assessing methylation changes of disease-specific genomic regions. With the mission to deliver the next level of simplicity, speed, and performance, Avida targeted methylation sequencing is based on a novel technology developed and perfected to preserve and retrieve methylation information from virtually any region of interest, even at very low DNA input levels. This chemistry is particularly optimized to detect circulating cell-free DNA (cfDNA), an emerging analyte that is proving increasingly important in molecular diagnostics. The technology also supports fragmented genomic DNA (gDNA) isolated from fresh/fresh-frozen samples or from FFPE-derived samples. With a fast turnaround and unparalleled sensitivity and specificity, Agilent's Avida targeted methylation sequencing technology can expertly advance your methylation-based liquid biopsy and epigenetics research.

Avida Target Enrichment Technology

At the core of the Avida target enrichment technology is an interlocked, three-dimensional structure, designed specifically for synergistic, indirect capture of intended DNA targets. In our proprietary design, a DNA scaffold forms with the intended target molecule when – and only when – more than one bridge probe is hybridized to the same target and stabilized by an anchor probe labeled with biotin (see [Figure 1](#)). Compared to traditional hybridization methods that use single and long biotinylated probes, the faster, more efficient Avida hybridization reaction is built on a much shorter, target-specific sequence used in the bridge probes. Consequently, a highly specific capture is achieved when the synergistic hybridization forms this interlocked structure and, in turn, binds to the streptavidin beads.

Avida probe panels are pre-formulated and ready-to-use to deliver consistent and best-in-class performance.

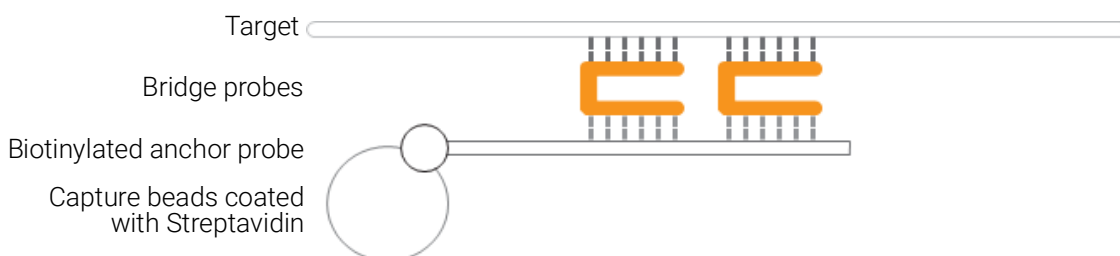


Figure 1 Hybridization of target to biotinylated anchor probe mediated by bridge probes

Overview of the Workflow

The starting material for the Avida Methyl workflow is cell-free DNA (cfDNA) or fragmented genomic DNA (gDNA) isolated from fresh/fresh-frozen samples or from FFPE-derived samples (refer to [“Input DNA samples”](#) on page 14 for detailed requirements on the DNA starting material).

The workflow features a PCR-free library preparation with a proprietary pre-amplification target capture process that maximizes capture efficiency, specificity, and uniformity. It also features our unique soft conversion – a next-generation bisulfite conversion method – to give you the most sample recovery possible. Altogether, this streamlined protocol delivers high conversion and library complexity which enables highly sensitive detection of minute methylation changes.

The workflow has four sections (see [Table 1](#)). You can expect the entire process to take approximately 6.5 hours to complete. The resulting libraries are ready for next-generation sequencing (NGS) on Illumina platforms.

Table 1 Workflow sections 1–4 and estimated time requirements for up to 8 samples

No.	Section	Description	Time
1	PCR-free library prep	The PCR-free library prep converts cfDNA or fragmented gDNA into adaptor-ligated molecules by end-repair and adaptor ligation	2 hours
2	Target capture	Target capture enriches specific targets from PCR-free libraries by probe hybridization and washes	1.5 hours
3	Soft conversion	Conversion includes DNA conversion, purification, and repair	1.5 hours
4	Indexing PCR and purification	Indexing PCR incorporates sample indexes and P5/P7 sequences necessary for Illumina sequencing	1.5 hours

Procedural and Safety Notes

Important practices for preventing contamination

The Avida Methyl technology is highly sensitive, making prevention of contamination of critical importance. Observe the following cautions when performing the workflow.

CAUTION

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.

CAUTION

To prevent PCR product contamination of samples throughout the workflow:

- 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. Reagents can be transferred from the pre-PCR area to the post-PCR area, but never from the post-PCR area to the pre-PCR area.
 - Maintain clean work areas. Before each assay, clean the surfaces that pose the highest risk of contamination using a solution of 10% bleach and water, or equivalent. Then, wipe the surfaces with water to remove bleach residue.
 - Always use dedicated pre-PCR pipettors with nuclease-free aerosol-resistant tips to pipette dedicated pre-PCR solutions.
 - Wear powder-free gloves. Use good laboratory hygiene, including changing gloves after contact with any potentially-contaminated surfaces.
-

CAUTION

To reduce the risk of contamination from an unconverted library to a converted library, change gloves at the specified steps in the protocol. Agilent strongly recommends separating the places where related equipment and supplies for conversion are located from the laboratory space where pre-PCR library prep and target capture is performed.

Safety notes

WARNING

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

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

Procedural notes

- This protocol has been developed using PCR 8-well strip tubes with individually attached caps or with tube cap strips. If using tube cap strips, 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.
- If handling multiple samples during washes (in [Chapter 4](#), “Target Capture”):
 - 1 Pre-aliquot reagents to a pipetting reservoir of appropriate size or to the tubes of a fresh strip tube.
 - 2 Use a multichannel pipette for washes.
- **Because the Avida Methyl workflow does not include a pre-capture amplification step, make sure no residual sample is left in the tube or pipette tip during sample transferring and pipetting to prevent any sample loss.**
- At certain steps that use magnetic beads, the complete removal of all supernatant is critical. The recommended technique for collecting residual supernatant at the bottom of the tubes depends on the magnet.
 - For magnets that collect beads on tube sides: Tap the magnet stand on the bench 5 times.
 - For magnets that collect beads at tube bottoms: Remove tubes from magnet and briefly spin. Return tubes to magnet before removing liquid.
- Reagents requiring -20°C storage are temperature sensitive and appropriate care should be taken during storage and handling. Avoid repeated freeze-thaw cycles. Make sure all reagents previously stored at -20°C are completely thawed before use.
- To process multiple samples, the protocol includes steps for preparation of reagent mixtures with overage, which are afterward distributed to the DNA library samples. To make reagent mixtures, scale reagents' volumes based on the number of reactions, using 10% excess volume to compensate for any pipetting loss. Add reagents in the order listed when preparing mixtures, then vortex to mix and briefly spin. Mixtures for preparation of 8 or 16 samples are shown in tables as examples.
- Possible stopping points, where samples may be stored at 4°C or -20°C , are marked in the protocol.

Materials Provided with Avida Methyl Kits and Panels

Table 2. lists the Avida Methyl reagent kits and the sub-kit boxes provided with each kit. The 16-reaction kits provide sufficient reagents for 2 assays with 8 samples. The 96-reaction kits provide sufficient reagents for 6 assays with 16 samples. The components of the sub-kit boxes are provided in **Chapter 8**, “Reagent Kit Contents,” starting on page 54.

Table 3 lists the Avida panel compatible with this Protocol.

Table 2 Avida Methyl reagent kits

Avida Methyl reagent kits	Sub-kit boxes	Storage Temperature
P/N G9419A Avida Methyl reagent kit with 1-16 index primer pairs for ILM, 16 reactions	Avida Methyl Reagent Box 1, 16 Reactions	–20°C
	Avida Methyl Reagent Box 2, 16 Reactions	Room temperature
	Avida Methyl and Duo Reagent Box 3, 16 Reactions	Room temperature
	Avida Beads Box, 16 Reactions	4°C
	Avida Index Primer Pairs 1–16 for ILM	–20°C
	Soft Conversion Reagent B*	Room temperature
P/N G9419B Avida Methyl reagent kit with 17-32 index primer pairs for ILM, 16 reactions	Avida Methyl Reagent Box 1, 16 Reactions	–20°C
	Avida Methyl Reagent Box 2, 16 Reactions	Room temperature
	Avida Methyl and Duo Reagent Box 3, 16 Reactions	Room temperature
	Avida Beads Box, 16 Reactions	4°C
	Avida Index Primer Pairs 17–32 for ILM	–20°C
	Soft Conversion Reagent B*	Room temperature
P/N G9429A Avida Methyl reagent kit with 1-96 index primer pairs for ILM, 96 reactions	Avida Methyl Reagent Box 1, 96 Reactions	–20°C
	Avida Methyl Reagent Box 2, 96 Reactions	Room temperature
	Avida Methyl and Duo Reagent Box 3, 96 Reactions	Room temperature
	Avida Beads Box, 96 Reactions	4°C
	Avida Index Primer Pairs 1–96 for ILM	–20°C
	Soft Conversion Reagent B*	Room temperature
P/N G9429B Avida Methyl reagent kit with 97-192 index primer pairs for ILM, 96 reactions	Avida Methyl Reagent Box 1, 96 Reactions	–20°C
	Avida Methyl Reagent Box 2, 96 Reactions	Room temperature
	Avida Methyl and Duo Reagent Box 3, 96 Reactions	Room temperature
	Avida Beads Box, 96 Reactions	4°C
	Avida Index Primer Pairs 97–192 for ILM	–20°C
	Soft Conversion Reagent B*	Room temperature

* Soft Conversion Reagent B may be shipped as part of the Avida Methyl and Duo Reagent Box 3 or it may be shipped in its own separate bag at room temperature.

Table 3 Compatible Avida Methyl panel

Capture Probes Panel	Target Size	Probes Size	Design ID	P/N for 16-reactions	P/N for 96-reactions
Avida Methyl 3400 DMR Cancer Panel	647 kb	743 kb	M3483261	5280-0058	5280-0059

NOTE

The **target size** is the size of the merged target regions used to design the capture probes for the panel. The **probes size** is the size of the merged genomic regions covered by the capture probes in the panel (see the *Covered.bed* file for the panel). You can download the bed files for a panel from the Agilent [SureDesign](#) website.

In the Avida Methyl 3400 DMR Cancer Panel, the differentially methylated regions used as the targets have a significant degree of overlap. Refer to the *Regions.bed* file for unmerged individual target annotation.

Online resource for Avida index sequences

Each member of the Avida Index Primer Pairs contains a unique 8-bp P5 or P7 index, resulting in dual-indexed NGS libraries. For complete index sequence information, download the Avida Index Sequence Resource Excel spreadsheet using this link: [Avida Index Resource](#).

Note that this link does not open a website. It automatically downloads the Excel spreadsheet to the default folder for downloaded files saved by your web browser. Locate the file in the folder to open it in Microsoft Excel or another compatible spreadsheet program. You can also find a link to the spreadsheet on the [Avida Methyl product page](#) on [www.agilent.com](#). The first tab of the spreadsheet provides instructions for use of the spreadsheet contents.

Additional Materials Required

Input DNA samples

The requirements for the DNA samples used as the starting material for the workflow are summarized in [Table 4](#). Input quantity ranges are based on Qubit Assay quantification of the DNA samples. DNA should be stored in low TE buffer (10 mM Tris-HCl, 0.1 mM EDTA, pH 8.0) or Qiagen Elution Buffer EB (10 mM Tris-HCl, pH 8.5). DNA needs to be at the appropriate size before you begin library preparation. For gDNA, Agilent recommends using Covaris shearing to achieve average insert sizes of 150–300 base pairs (see [Chapter 2](#), “gDNA Fragmentation” for instructions). Since cfDNA typically has an average size of 170 base pairs, no further fragmentation is required.

Table 4 Starting material requirements by DNA sample type

Sample type	Input quantity	Fragmentation requirements	Storage buffer
cfDNA	3–100 ng [*]	No fragmentation required as average fragment size is already 170 bp	Low TE buffer or Qiagen Buffer EB
gDNA from fresh or fresh-frozen sample or FFPE-derived sample	10–100 ng [†]	Fragmentation with Covaris shearing to insert sizes of 150–300 bp	Low TE buffer or Qiagen Buffer EB

^{*} Libraries can be generated with inputs of cfDNA as low as 3 ng. For detection of methylation changes, increasing the input improves performance.

[†] For FFPE-derived gDNA, use the maximum amount of sample available in this range.

NOTE

The Avida workflow is not compatible with the Seraseq ctDNA Complete Reference Material or the Seraseq ctDNA Reference Material V2 available from SeraCare. Contact Agilent [Technical Support](#) with questions.

Additional reagents and equipment required

See [Table 5](#) through [Table 7](#) for additional reagents and equipment required to complete the workflow.

Table 5 Additional required reagents

Description	Vendor and Part Number
100% Ethanol (Ethyl Alcohol, 200 proof)	Millipore p/n EX0276, or equivalent
1X Low TE Buffer (10 mM Tris-HCl, pH 8.0, 0.1 mM EDTA)	Thermo Fisher Scientific p/n 12090-015, or equivalent
Qubit dsDNA HS or BR Assay Kit (optional) 100 assays 500 assays	Thermo Fisher Scientific p/n Q32851 (HS) or Q32850 (BR) p/n Q32854 (HS) or Q32853 (BR)
Purification Beads: AMPure XP for PCR purification 5 mL 60 mL	Beckman Coulter Genomics p/n A63880 p/n A63881

Table 6 Required equipment

Description	Vendor and Part Number
Thermal Cycler with 96-well, 0.2 mL block	Various suppliers
PCR 8-well strip tubes, 0.2-mL, with individually attached caps or with strip caps	VWR p/n 76318-798, Genesee p/n 27-125U, or consult the thermal cycler manufacturer's recommendations
Nucleic acid analysis system (instrument and consumables)	Select one system from Table 7 on page 15
Qubit 4 Fluorometer and Qubit Assay Tubes (optional)	Thermo Fisher Scientific p/n Q33238 and p/n Q32856
Heat blocks or water baths heated to 37°C and 50°C	General laboratory supplier
DNA LoBind Tubes, 1.5-mL PCR clean, 250 pieces	Eppendorf p/n 022431021 or equivalent
Microcentrifuge	Eppendorf microcentrifuge, model 5417C or equivalent
Centrifuge for 8-well strip tubes, 3000–5000 rpm	General laboratory supplier
Centrifuge for 96-well plate (if using 96-reaction kits)	General laboratory supplier
DynaMag-96 Side Magnet or Magnum EX Universal 96-well Magnet Plate	Thermo Fisher Scientific p/n 12331D or Alpaqua p/n A000380
Magnet for 1.5-mL tubes (optional)	Thermo Fisher Scientific p/n 12321D or equivalent
Multichannel and single channel pipettes	Rainin Pipet-Lite Multi Pipette or equivalent
Reagent reservoirs for use with multichannel pipettes (optional)	VWR 10015-236 or equivalent
Sterile, nuclease-free aerosol barrier pipette tips, vortex mixer, cold rack, ice bucket, and powder-free gloves	General laboratory supplier

Table 7 Nucleic acid analysis platform options – select one

Analysis System	Vendor and Part Number Information
Agilent 4200/4150 TapeStation Instrument	Agilent p/n G2991AA/G2992AA
Consumables:	
96-well sample plates	p/n 5042-8502
96-well plate foil seals	p/n 5067-5154
8-well tube strips	p/n 401428
8-well tube strip caps	p/n 401425
D1000 ScreenTape	p/n 5067-5582
D1000 Reagents	p/n 5067-5583
Agilent 5200/5300/5400 Fragment Analyzer Instrument	Agilent p/n M5310AA/M5311AA/M5312AA
Consumables:	
<i>Select from the following kits*</i>	
NGS Fragment Kit (1-6000 bp)	p/n DNF-473-0500
Small Fragment Kit (1-1500 bp)	p/n DNF-476-0500
HS NGS Fragment Kit (1-6000 bp)	p/n DNF-474-0500
HS Small Fragment Kit (1-1500 bp)	p/n DNF-477-0500

* Consider factors such as sample type and quality, expected target insert size, and sample input quantity when selecting a DNA Analysis Kit for the Fragment Analyzer system.

Equipment required for gDNA fragmentation

[Table 8](#) lists information on equipment for fragmentation of gDNA samples using mechanical shearing. Sample fragmentation is required for high-quality gDNA samples and FFPE-derived gDNA samples but not for cfDNA samples.

Table 8 Equipment for mechanical shearing of gDNA samples

Description	Vendor and Part Number	Usage Notes
Covaris Sample Preparation System	Covaris model E220	Additional Covaris instrument models and sample holders may be used after optimization of shearing conditions.
Covaris microTUBE sample holders	Covaris p/n 520045	

2

gDNA Fragmentation

Step 1. Prepare for shearing [18](#)

Step 2. Shear the gDNA samples [18](#)

This chapter describes the steps to fragment gDNA samples using mechanical shearing. gDNA samples must be fragmented prior to starting the Avida Methyl workflow.

NOTE

Fragmentation is only required for gDNA samples. If you are using cfDNA samples, proceed to [Chapter 3](#), "PCR-Free Library Preparation."

Step 1. Prepare for shearing

NOTE

The shearing instructions provided here were optimized using a Covaris model E220 instrument and the 130- μ L Covaris microTUBE. Consult the manufacturer's recommendations for use of other Covaris instruments or sample holders to achieve the same target gDNA fragment size of 150–300 bp.

- 1 Set up the Covaris instrument according to the manufacturer's instructions. Allow enough time (typically 30–60 minutes) for instrument degassing and water bath chilling before starting the protocol.
- 2 Dilute 10–100 ng of each gDNA sample with 1X low TE buffer to a final volume of 50 μ L. Vortex well to mix, then briefly spin. Keep the samples on ice.

Step 2. Shear the gDNA samples

- 1 For each gDNA sample, transfer the 50- μ L sample into a Covaris microTUBE.
- 2 Spin the microTUBE for 30 seconds to collect the liquid and to remove any bubbles from the bottom of the tube.
- 3 Secure the microTUBE in the tube holder and shear the gDNA with the settings in [Table 9](#).

Table 9 Shear settings for Covaris E-series instrument (SonoLab software v7 or later)

Setting	High-quality gDNA	FFPE-derived gDNA
Duty Factor	10%	10%
Peak Incident Power (PIP)	175	175
Cycler per Burst	200	200
Treatment Time	2 \times 60 seconds (see two-round instructions below)	240 seconds
Bath Temperature	2° to 8°C	2° to 8°C

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

- a Shear for 60 seconds
 - b Spin the microTUBE for 10 seconds
 - c Vortex the microTUBE at high speed for 5 seconds
 - d Spin the microTUBE for 10 seconds
 - e Repeat [step a](#) through [step d](#)
- 4 After shearing, put the Covaris microTUBE back into the loading and unloading station.
 - 5 Spin the microTUBE briefly. While keeping the snap-cap on, insert a pipette tip through the pre-split septum, then slowly remove the sheared gDNA.
 - 6 Transfer the sheared gDNA sample (approximately 50 μ L) to a strip tube sample well. Keep the samples on ice.

2 gDNA Fragmentation

Step 2. Shear the gDNA samples

- 7 After transferring the gDNA sample, spin the microTUBE briefly to collect any residual sample volume. Transfer any additional collected liquid to the sample well used in [step 6](#) above.

The 50-μL sheared gDNA samples are now ready for library preparation. Proceed directly to [Chapter 3](#), “PCR-Free Library Preparation.”

This is not a stopping point.

3

PCR-Free Library Preparation

Step 1. End prep (end repair and dA-tailing) [21](#)

Step 2. Ligation [23](#)

Step 3. Library bead binding [25](#)

This chapter describes the steps for the PCR-free library preparation section of the Avida Methyl workflow. NGS libraries are prepared using adaptors that include Unique Molecular Identifiers (UMIs).

See [Table 4](#) on page 14 for DNA sample input requirements.

This chapter uses the reagents listed in [Table 10](#).

Table 10 Reagents for PCR-free library preparation

Reagent	Usage Notes	Kit/Storage Location
End Prep Buffer (tube with purple cap)	Thaw at room temperature; vortex before use.	Avida Methyl Reagent Box 1, stored at –20°C
End Prep Enzyme (tube with blue cap)	Place on ice or cold block just before use.	
Ligation Buffer (tube with green cap or bottle)	Thaw at room temperature; vortex before use.	
Ligation Enzyme (tube with yellow cap)	Place on ice or cold block just before use.	
Adaptor for ILM (tube with orange cap)	Thaw at room temperature then keep on ice.	
Hyb Blocker (tube with red cap)	Thaw at room temperature.	
Library Binding Beads (tube with white cap or bottle)	Equilibrate to room temperature for at least 15 minutes.	Avida Beads Box, stored at 4°C
Nuclease-free Water (tube with clear cap or bottle)	—	Avida Methyl Reagent Box 2, stored at Room Temperature
Library Wash Buffer (bottle)	—	
Avida Methyl 3400 DMR Cancer Panel	Thaw at room temperature then keep on ice.	Stored at –20°C

CAUTION

Review the “[Procedural and Safety Notes](#)” on page 10 before getting started.

Because the Avida Methyl workflow does not include a pre-capture amplification step, make sure no residual sample is left in the tube or pipette tip during sample transferring and pipetting to prevent any sample loss.

NOTE

Prior to starting a workflow, begin preheating the stock bottle of Hyb Wash Buffer 1 to ensure that it is sufficiently heated and solubilized prior to its use in [Chapter 4](#). The Hyb Wash Buffer 1 is provided in the Avida Methyl Reagent Box 2.

- If a water bath is available, heat the stock bottle in a 50°C water bath for 10 minutes. Then, transfer the bottle to a heat block set to 50°C. Leave the bottle on the 50° heat block until the buffer is used in [Chapter 4](#).
- If a water bath is not available, place the stock bottle on a heat block set to 50°C. Leave the bottle on the 50°C heat block until the buffer is used in [Chapter 4](#).

If the bottle does not fit in the heat block, place the bottle on top of the heat block, setting it on its side to increase surface area contact. A thermal cycler thermal block set to 50°C may be used in place of a heat block.

Step 1. End prep (end repair and dA-tailing)

- 1 Thaw the End Prep Buffer at room temperature, then thoroughly vortex. Inspect the buffer to make sure that no small particles are visible. If particles are visible, continue to vortex until they dissolve.
- 2 For each sample, add an appropriate quantity of DNA (based on sample type, see below) to a strip tube in a volume of 50 µL. If the sample volume is <50 µL, add Nuclease-free Water to bring the volume to 50 µL. See [Table 4](#) on page 14 for complete information on the DNA sample input requirements. Keep DNA samples on ice.
 - cfDNA samples: 3–100 ng
 - fragmented gDNA samples: 10–100 ng (prepared in [Chapter 2](#))
- 3 Prepare the appropriate volume of End Prep Mix by combining the reagents in [Table 11](#) at room temperature. Combine the reagents in the order shown. Gently vortex the End Prep Mix at low speed, then briefly spin.

Once prepared, keep the End Prep Mix on ice or use it immediately. The End Prep Mix is stable for up to 2 hours at 4°C.

Table 11 Preparation of End Prep Mix

Reagent	Volume for 1 reaction	Volume for 8 reactions (includes excess)	Volume for 16 reactions (includes excess)
End Prep Buffer	7 µL	63 µL	126 µL
End Prep Enzyme	3 µL	27 µL	54 µL
Total	10 µL	90 µL	180 µL

- 4 Add 10 µL of End Prep Mix to each tube containing a DNA sample. Mix by gentle vortexing or by pipetting up and down 15–20 times using a pipette set to 50 µL. Briefly spin the strip tubes then keep on ice.

A small amount of bubbles in the end prep reactions does not affect performance.

PCR-Free Library Preparation**Step 1. End prep (end repair and dA-tailing)**

- 5 Program the thermal cycler to run the end prep thermal cycler program shown in [Table 12](#). Set the heated lid to 75°C. Load the strip tubes and run the program.

Table 12 Thermal cycler program for end prep

Step	Temperature	Time
Step 1	20°C	30 minutes
Step 2	65°C	30 minutes
Step 3	4°C	Hold

- 6 Start thawing the Ligation Buffer and the Adaptor for ILM at room temperature in preparation for the next step. Thoroughly vortex once thawed. Keep the thawed Adaptor for ILM on ice. Inspect the Ligation Buffer to make sure that no small particles are visible. If particles are visible, continue to vortex until they dissolve.

Step 2. Ligation

- 1 While the end prep thermal cycler program is running, prepare the Ligation Master Mix by combining the reagents in [Table 13](#) in a 1.5-mL tube at room temperature. Vortex the Ligation Master Mix, then briefly spin the tube. Once prepared, keep it on ice until needed or use it immediately.

CAUTION

The Ligation Buffer and Ligation Enzyme are viscous. Take care when pipetting these reagents and mixtures containing these reagents.

The Ligation Master Mix is stable for up to 2 hours at 4°C.

Table 13 Preparation of Ligation Master Mix

Reagent	Volume for 1 reaction	Volume for 8 reactions (includes excess)	Volume for 16 reactions (includes excess)
Ligation Buffer	25 µL	225 µL	450 µL
Ligation Enzyme	6 µL	54 µL	108 µL
Total	31 µL	279 µL	558 µL

- 2 When the end prep thermal cycler program is complete, remove the strip tubes containing the end-repaired/dA-tailed DNA samples. Keep at room temperature.

Make sure to stop the thermal cycler program to allow the thermal block and heated lid to equilibrate to room temperature.

- 3 Add 5 µL of Adaptor for ILM to each DNA sample.

CAUTION

Do NOT add Adaptor for ILM to the Ligation Master Mix. Avoid exposing the Adaptor to room temperature conditions.

- 4 Add 31 µL of Ligation Master Mix to each sample. Mix by gentle vortexing or by pipetting up and down 15–20 times using a pipette set to 70 µL. Briefly spin the samples.
- 5 Make sure that the thermal cycler has equilibrated to room temperature. Program the thermal cycler to run the ligation thermal cycler program shown in [Table 14](#) with the heated lid off. Load the strip tubes and run the program.

Table 14 Thermal cycler program for ligation

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

- 6 During the thermal cycler program for ligation, prepare the following reagents as described below in preparation for upcoming steps.
- Equilibrate the Library Binding Beads to room temperature for at least 15 minutes before use in [“Step 3. Library bead binding.”](#)
 - Heat the Hyb Buffer to 37°C in a water bath or heat block for at least 10–20 minutes, or until the Hyb Buffer is completely solubilized. If the Hyb Buffer is supplied in a bottle, and the bottle does not fit into the heat block, place the bottle on its side (to increase surface area contact) on top of the heat block. Once the Hyb Buffer is completely solubilized, transfer to room temperature until use. The Hyb Buffer is used in [Chapter 4](#), “Target Capture.”

NOTE

If your laboratory does not have a water bath or heat block, you can heat the tube or bottle of Hyb Buffer by placing it on top of the thermal block of a thermal cycler set to 37°C. More time may be required to completely solubilize the buffer with this technique, especially if the Hyb Buffer is supplied in a bottle.

- Thaw the Hyb Blocker, Hyb Enhancer, and Avida Panel at room temperature. Transfer the Avida Panel to ice once thawed. Keep the Hyb Blocker and Hyb Enhancer at room temperature until use. The Hyb Blocker and Avida Panel are used to prepare Hyb Mix 1 in [Table 15](#). The Hyb Enhancer is used in [Chapter 4](#), “Target Capture.”

CAUTION

The Hyb Enhancer must be thawed at room temperature. Once thawed, thoroughly vortex the Hyb Enhancer to solubilize any precipitates.

Step 3. Library bead binding

- 1 Once the thermal cycler program for ligation completes, remove the strip tubes containing the adaptor-ligated DNA samples from the thermal cycler.
- 2 Thoroughly vortex the stock bottle or stock tube of Library Binding Beads to mix.

Note that some bead aggregates may remain even after vortexing. The aggregates do not impact performance.
- 3 Add 87 µL (~0.9X volume) of Library Binding Beads to each adaptor-ligated DNA sample, bringing the sample volume to 183 µL. Mix by pipetting up and down 15–20 times using a pipette set to 150 µL. Spin for 1 to 2 seconds to collect the liquid at the bottom of the tube, making sure the beads do not settle.

Pipette slowly to avoid creating excess bubbles. Bubbles can compromise library binding efficiency.
- 4 Incubate the library bead binding reactions for 10 minutes at room temperature.
- 5 **During the 10-minute incubation**, prepare the Hyb Mix 1 by combining the reagents in [Table 15](#) in a 1.5-mL tube. Vortex the Hyb Mix 1 thoroughly, then briefly spin the tube. Set aside at room temperature until needed in the next chapter.

NOTE

When handling multiple samples, Agilent recommends aliquoting 26 µL of Hyb Mix 1 into the tubes of a fresh strip tube (one tube per sample). Then, during [step 1](#) on [page 28](#), use a multichannel pipette to transfer 24 µL of Hyb Mix 1 from the strip tube to the beads bound with adaptor-ligated library. This technique helps prevent the beads from drying out.

Table 15 Preparation of Hyb Mix 1

Reagent	Volume for 1 reaction	Volume for 8 reactions (includes excess)	Volume for 16 reactions (includes excess)
Nuclease-Free Water	17.5 µL	157.5 µL	315 µL
Avida Methyl Panel	4 µL	36 µL	72 µL
Hyb Blocker	2.5 µL	22.5 µL	45 µL
Total	24 µL	216 µL	432 µL

- 6 When the 10-minute incubation is complete, place the strip tubes on a magnet for at least 2 minutes or until the solution is clear.

NOTE

The solution may take longer than 2 minutes to become clear. Visually inspect the tubes to make sure that the beads have fully aggregated and the solution is clear before proceeding.

- 7 Remove and discard all supernatant from each tube without removing or disturbing the beads. **Leave the strip tubes on the magnet until instructed to remove them.**

PCR-Free Library Preparation**Step 3. Library bead binding**

- 8** Add 180 μ L of Library Wash Buffer to each tube without disturbing the bead pellet. Incubate for 2 minutes at room temperature.
- 9** Remove and discard all supernatant from each tube.
- 10** (Optional) Repeat [step 8](#) through [step 9](#) for an additional wash.

This second wash may help improve percent on-target in sequencing results, but it may slightly reduce molecular recovery.

- 11** Collect all residual supernatant at the bottom of the tubes, then remove and discard the residual supernatant from each tube using a 20 μ L pipette.

Refer to [“Procedural notes”](#) on page 11 for best practices on removing all traces of supernatant based on magnet type.

- 12** Remove the strip tubes from the magnet.

The adaptor-ligated libraries remain bound to the Library Binding Beads.

- 13** Proceed **immediately** to [step 1](#) on [page 28](#) (in [Chapter 4](#), “Target Capture”) and add the Hyb Mix 1 that you prepared earlier to make sure that the beads do not dry out.

4 Target Capture

- Step 1. Hybridization [28](#)
 Step 2. Capture Bead Preparation [29](#)
 Step 3. Bead Capture [31](#)
 Step 4. Washing (Bead-Bound DNA Hybrids) [32](#)

This chapter describes the steps for the target capture section of the Avida Methyl workflow. The adaptor-ligated libraries are hybridized with a target-specific panel. After hybridization, the targeted molecules are captured on Capture Beads coated with Streptavidin.

This chapter uses the reagents listed in [Table 16](#).

Table 16 Reagents for target capture

Reagent	Usage Notes	Kit/Storage Location
Hyb Buffer (tube with clear cap or bottle)	Heat to 37°C in water bath or heat block for at least 10–20 minutes before use. Once solution becomes clear, keep at room temperature.	Avida Methyl Reagent Box 1, stored at –20°C
Hyb Enhancer (amber tube with green cap)	Thaw at room temperature.	
Capture Beads (tube with amber cap)	Thoroughly vortex to mix before use.	Avida Beads Box, stored at 4°C
Nuclease-free Water (tube with clear cap or bottle)	—	Avida Methyl Reagent Box 2, stored at Room Temperature
Hyb Wash Buffer 1 (bottle)	Heat the stock bottle at 50°C in water bath or heat block until use.	
Hyb Wash Buffer 2 (bottle)	—	
Hyb Mix 1	Prepared in previous chapter (page 25) and kept at room temperature until use.	

Step 1. Hybridization

- 1 Add 24 μL of Hyb Mix 1 (prepared on [page 25](#)) to the beads bound with adaptor-ligated library generated in the last step of the previous chapter. Resuspend the beads by gentle vortexing or by pipetting up and down 15–20 times. Briefly spin the strip tubes.

NOTE

If during preparation of Hyb Mix 1 ([page 25](#)) you distributed 26 μL of Hyb Mix 1 into the tubes of a fresh strip tube, you can now use a multichannel pipette to transfer 24 μL from the strip tube to the beads bound with adaptor-ligated library.

- 2 Thoroughly vortex the Hyb Buffer that has been kept at room temperature.
If a precipitate is visible in the Hyb Buffer, reheat to 37°C (as described on [page 24](#)) to resolubilize prior to use.
- 3 Prepare the Hyb Mix 2 by combining the reagents in [Table 17](#) in a 1.5-mL tube. Vortex the Hyb Mix 2 thoroughly, then briefly spin the tube. Keep at room temperature.

Keep the Hyb Buffer at room temperature until later use.

Table 17 Preparation of Hyb Mix 2

Reagent	Volume for 1 reaction	Volume for 8 reactions (includes excess)	Volume for 16 reactions (includes excess)
Hyb Buffer	30 μL	270 μL	540 μL
Hyb Enhancer	6 μL	54 μL	108 μL
Total	36 μL	324 μL	648 μL

- 4 Add 36 μL of Hyb Mix 2 to each bead mix containing Hyb Mix 1 (from [step 1](#) above), bringing the total volume of each hybridization reaction to 60 μL .
- 5 Mix by pipetting up and down 15–20 times using a pipette set to 50 μL . Spin for 1 to 2 seconds to collect the liquid at the bottom of the tube, making sure the beads do not settle.
- 6 Program the thermal cycler to run the hybridization thermal cycler program shown in [Table 18](#). Set the heated lid to 103°C. Load the strip tubes and run the program.

Table 18 Thermal cycler program for hybridization

Step	Temperature	Time
Step 1	98°C	2 minutes
Step 2	$\geq 2.5^\circ\text{C}/\text{second}$ ramp down to 60°C	
Step 3	60°C	60 minutes
Step 4	60°C	Hold

NOTE

The thermal cycler program in [Table 18](#) uses a 60-minute hybridization which is sufficient for most applications. A longer hybridization time (up to 16 hours) may improve capture efficiency, especially for panels >500 kb.

Step 2. Capture Bead Preparation

CAUTION

Hyb Wash Buffer 1 may become cloudy or form crystals during storage. Make sure that the solution solubilized during the 50°C heating. Thoroughly vortex or invert the stock bottle before use.

Capture bead preparation can be performed during the last 15 minutes of the hybridization thermal cycler program.

- 1 Thoroughly vortex the Hyb Buffer that has been kept at room temperature since its last use in [“Step 1. Hybridization.”](#)
- 2 Thoroughly pipette or vortex the Capture Beads stock to make sure the solution is homogeneous.
- 3 8 µL of Capture Beads are required per reaction. Calculate the total volume of beads needed for all the samples, including 10% overage, and transfer that volume into a 1.5-mL tube.

NOTE

Preparing the Capture Beads in a 1.5-mL tube requires a magnet that accommodates that tube size. If needed, the required volume of Capture Beads can be divided among multiple PCR tubes.

- 4 Place the tube of Capture Beads on the magnet for 1 minute or until the solution becomes clear. Remove and discard the supernatant. Remove the tube from the magnet.
- 5 Add the appropriate volume of the preheated Hyb Wash Buffer 1 to the Capture Beads using the guidelines in [Table 19](#). Mix well by vortexing or by pipetting up and down at least 15–20 times.

Make sure the beads are fully resuspended and well mixed. When working with wash volumes >100 µL, mixing by pipetting is more effective than vortexing.

Table 19 Wash volumes of Hyb Wash Buffer 1 based on number of hybridization reactions

Number of reactions	Volume of Hyb Wash Buffer 1
1–16	100 µL
17–32	200 µL
33–48	300 µL
49–64	400 µL
65–80	500 µL
81–96	600 µL

- 6 Briefly spin the tube of Capture Beads, then place the tube on the magnet for 1 minute or until the solution becomes clear. Remove and discard the supernatant.
- 7 Perform an additional wash by repeating [step 5](#) through [step 6](#) for a total of two washes.
Place the Hyb Wash Buffer 1 back on the 50°C heat block or water bath.

Target Capture**Step 2. Capture Bead Preparation**

- 8 Remove the tube of Capture Beads from the magnet. Resuspend the beads in Hyb Buffer at the same volume you originally started with (as calculated in [step 3](#), above). Mix by pipetting up and down 15–20 times (avoid creating excess bubbles).

NOTE

Make sure to use Hyb Buffer to resuspend the Capture Beads.

-
- 9 Add 8 μ L of resuspended Capture Beads to each tube of a fresh strip tube (one for each hybridization reaction).

Step 3. Bead Capture

- 1 Once the hybridization thermal cycler program completes, remove the strip tubes containing the hybridization reactions from the thermal cycler and set them aside. Stop the hybridization thermal cycler program, and start the washing thermal cycler program shown in [Table 20](#). Set the heated lid to 75°C.

Table 20 Thermal cycler program for washing

Step	Temperature	Time
Step 1	60°C	Hold

- 2 Briefly spin the strip tubes then place on a magnet. Keep the tubes on the magnet for 30 seconds or until the solution becomes clear (up to 1 minute).
- 3 **Carefully transfer the supernatant from each tube** (making sure to get all the supernatant) to a tube containing an 8-μL aliquot of prepared Capture Beads (from [step 9](#) on [page 30](#)). Mix by pipetting up and down 15–20 times (avoid vortexing at this step). Spin for 1 to 2 seconds to collect the liquid at the bottom of the tubes, making sure the beads do not settle.
- 4 Perform bead binding by placing the strip tubes on the thermal cycler running at 60°C for 10 minutes.
- 5 After 10 minutes, remove the strip tubes from the thermal cycler and place on a magnet for 30 seconds or until the solution becomes clear (up to 1 minute). Remove and discard all supernatant.

Leave the thermal cycler running at 60°C.
- 6 Remove the strip tubes from the magnet and immediately proceed to [“Step 4. Washing \(Bead-Bound DNA Hybrids\).”](#)

Step 4. Washing (Bead-Bound DNA Hybrids)

When handling multiple samples, Agilent recommends transferring Hyb Wash Buffer 1 and Hyb Wash Buffer 2 into reagent reservoirs and pipetting the buffers using a multichannel pipette.

Refer to “[Procedural notes](#)” on page 11 for best practices on removing all traces of supernatant based on magnet type.

CAUTION

For [step 1](#) through [step 5](#) in this section, do NOT vortex or spin the samples. Carefully handle the tubes to prevent any splashing.

- 1 Add 150 μ L of preheated Hyb Wash Buffer 1 to each tube containing the Capture Beads/library DNA hybrids. Mix by pipetting up and down at least 15–20 times, making sure the beads are fully resuspended.
- 2 Put the strip tubes on a magnet for 30 seconds or until the solution becomes clear. Remove and discard the supernatant from each tube. Use a 20 μ L pipette to remove all residual Hyb Wash Buffer 1.
- 3 Remove the strip tubes from the magnet. Add 100 μ L of Hyb Wash Buffer 1 to each tube. Mix by pipetting up and down 15–20 times, making sure the beads are fully resuspended. *(Complete resuspension is critical.)* **Then, transfer the entire mix (beads, buffer, and any bubbles that may have formed) to a fresh strip tube.**
- 4 Place the strip tubes on the thermal cycler running at 60°C for 3 minutes.
- 5 After 3 minutes, remove the strip tubes from the thermal cycler and stop the thermal cycler wash program.
- 6 Place the strip tubes on the magnet for 30 seconds or until the solution becomes clear. Remove and discard the supernatant from each tube. Use a 20 μ L pipette to remove all residual Hyb Wash Buffer 1.

Promptly execute the next steps ([step 7](#) and [step 8](#)) to avoid over-drying the bead pellets.
- 7 Remove the strip tubes from the magnet.
- 8 Add 150 μ L of room temperature Hyb Wash Buffer 2 to each tube. Mix by pipetting up and down at least 15–20 times, making sure the beads are fully resuspended. *(Complete resuspension is critical.)* Briefly spin. Proceed to [Chapter 5](#), “Soft Conversion and Repair.”

Stopping Point If you do not continue to the next step, store the samples in Hyb Wash Buffer 2 at 4°C overnight.

5 Soft Conversion and Repair

- Step 1. Soft Conversion [34](#)
 Step 2. Bead Purification [36](#)
 Step 3. Washing (Bead-Bound Converted Libraries) [37](#)
 Step 4. DNA Repair [38](#)

This chapter describes the steps for the soft conversion and repair section of the Avida Methyl workflow. The adaptor-ligated libraries are treated with soft conversion reagents to deaminate any unmethylated cytosines, thereby converting them to uracils. The converted DNA is then purified, washed, and repaired. During the PCR indexing step ([Chapter 6](#), "Indexing PCR, Library Purification, and Quality Assessment") the uracils are replaced with thymines in the PCR products.

This chapter uses the reagents listed in [Table 21](#).

Table 21 Reagents for soft conversion and repair

Reagent	Usage Notes	Kit/Storage Location
Soft Conversion Reagent A (amber tube with amber cap)	Thaw at room temperature; vortex to mix.	Avida Methyl Reagent box 1; stored at –20°C
Soft Conversion Reagent B (amber tube with blue cap)	Vortex to mix; discard tube after use (do not re-use).	Stored at room temperature; either included with Avida Methyl and Duo Reagent Box 3 or shipped separately
Soft Conversion Binding Beads (bottle)	Vortex to mix.	Avida Methyl and Duo Reagent Box 3, stored at room temperature
5X Soft Conversion Wash Buffer (bottle)	Add ethanol prior to initial use.	
Soft Conversion Elution Buffer (tube with green cap)	—	
Soft Conversion Repair Solution (tube with yellow cap)	—	
Nuclease-Free Water (tube with clear cap or bottle)	—	
100% ethanol	—	Not provided

Step 1. Soft Conversion

CAUTION

Agilent recommends designating a separate area for conversion work to avoid contamination with unconverted materials. Always change to a new pair of gloves for conversion work.

- 1 Briefly spin the strip tubes containing the Capture Beads/library DNA hybrids in Hyb Wash Buffer 2 (from the last step of [Chapter 4](#)).
- 2 Place the strip tubes on the magnet for 1 minute or until the solution becomes clear. Remove and discard the supernatant from each tube.
- 3 Collect all residual supernatant at the bottom of the tubes, then remove and discard the residual supernatant from each tube using a 20 μ L pipette. Immediately remove the tubes from the magnet.

CAUTION

Removing all residual Hyb Wash Buffer 2 from the beads is critical. Residual buffer may lower the conversion rate.

Do not leave the tubes on the magnet once the buffer is completely removed. Leaving dry beads on the magnet could cause the beads to aggregate, resulting in low recovery yield.

- 4 Add 11.5 μ L of Soft Conversion Elution Buffer to each tube. Gently pulse vortex to ensure the beads are fully resuspended in the solution. Spin the strip tubes for 1–2 seconds to collect the liquid at the bottom of the tube, making sure the beads do not settle.

NOTE

For gentle pulse vortexing, reduce the vortex speed to 50% power, then pulse vortex the tubes for 1–2 seconds at a time for a total of 10 times.

NOTE

When handling multiple samples, Agilent recommends aliquoting 12.5 μ L of Soft Conversion Elution Buffer into the tubes of a fresh strip tube (one tube per sample). Then, use a multichannel pipette to transfer 11.5 μ L of the buffer from the strip tube to the beads bound with adaptor-ligated library. This technique helps prevent the beads from drying out.

- 5 Incubate the tubes at room temperature for 10 minutes.
- 6 During the 10-minute incubation, prepare the Soft Conversion Mix by combining the components in [Table 22](#) in a 1.5-mL tube.

Thoroughly vortex Soft Conversion Reagent A and Soft Conversion Reagent B before adding them to the mix. It is normal to see some white crystals forming around the tube and lid of Soft Conversion Reagent B.

NOTE

Each tube of Soft Conversion Reagent B is provided as a single-use reagent. Once the tube is opened, the solution is susceptible to oxidization. Discard the tube after use.

Soft Conversion and Repair

Step 1. Soft Conversion

Vortex the Soft Conversion Mix thoroughly, then briefly spin the tube. The Soft Conversion Mix needs to be prepared fresh each time and used within 10 minutes after preparation. Keep the cap of the tube closed as much as possible.

Table 22 Preparation of Soft Conversion Mix

Reagent	Volume for 1 reaction	Volume for 8 reactions (includes excess)	Volume for 16 reactions (includes excess)
Soft Conversion Reagent A	3.5 μ L	31.5 μ L	63 μ L
Soft Conversion Reagent B	20 μ L	180 μ L	360 μ L
Total	23.5 μL	211.5 μL	423 μL

- 7 After the 10-minute incubation ([step 5](#), above), place the strip tubes on the magnet for 1 minute or until the solution becomes clear (up to 2 minutes).
- 8 **Carefully transfer the supernatant from each tube of the strip tube** (making sure to get all the supernatant) to a fresh strip tube.
- 9 Add 23.5 μ L of freshly prepared Soft Conversion Mix to each tube. Mix by pipetting up and down 15–20 times then immediately close the cap.
- 10 Program the thermal cycler to run the conversion thermal cycler program shown in [Table 23](#). Set the heated lid to 85°C. Load the strip tubes and run the program.

Table 23 Thermal cycler program for conversion

Step	Temperature	Time
Step 1	75°C	25 minutes *
Step 2	10°C	Hold

* If conversion >99.5% is desired for your assay, increase the duration of Step 1 from 25 minutes to 35 minutes. While this added duration maximizes conversion, it will reduce molecular recovery by 10–15%.

CAUTION

As soon as the thermal cycler program reaches the 10°C hold step, proceed immediately to [“Step 2. Bead Purification.”](#) Prolonged holding at 10°C may lower the recovery yield.

Step 2. Bead Purification

- 1 Once the thermal cycler program for conversion reaches the 10°C hold step, remove the strip tubes from the thermal cycler.

NOTE

You may see a brown coating in the tubes following conversion. This appearance is normal.

- 2 Thoroughly vortex the stock of Soft Conversion Binding Beads, then add 150 µL of beads to each of the tubes containing a converted library. Mix by pipetting up and down 15–20 times, making sure the beads are fully resuspended. Spin for 1 to 2 seconds to collect the liquid at the bottom of the tubes, making sure the beads do not settle.
- 3 Incubate the tubes at room temperature for 10 minutes.
- 4 During the 10-minute incubation, prepare a 0.1X dilution of the Soft Conversion Elution Buffer by combining the components in [Table 24](#) in a 1.5-mL tube.

The 0.1X Soft Conversion Elution Buffer is used in [“Step 4. DNA Repair”](#) on page 38.

Table 24 Preparation of 0.1X Soft Conversion Elution Buffer

Reagent	Volume for 1 reaction	Volume for 8 reactions (includes excess)	Volume for 16 reactions (includes excess)
Soft Conversion Elution Buffer	1.5 µL	16.5 µL	33 µL
Nuclease-Free Water	13.5 µL	148.5 µL	297 µL
Total	15 µL	165 µL	330 µL

- 5 After the 10-minute incubation ([step 3](#), above), carefully place the strip tubes on the magnet for 30 seconds or until the solution becomes clear (up to 1 minute). Use a 20 µL pipette to remove and discard all supernatant.
- 6 Remove the tubes from the magnet and proceed immediately to [“Step 3. Washing \(Bead-Bound Converted Libraries\).”](#)

Step 3. Washing (Bead-Bound Converted Libraries)

NOTE

Before using the Soft Conversion Wash Buffer for the first time, prepare a 1X stock by adding 100% ethanol to the stock bottle containing 5X Soft Conversion Wash Buffer. Vortex to mix, then mark the check box on the bottle to indicate that ethanol was added. Keep bottle tightly capped to avoid ethanol evaporation.

- The bottle of 5X Soft Conversion Wash Buffer supplied with the 16-reaction kits (p/n 5271-0115) requires 8 mL of 100% ethanol.
- The bottle of 5X Soft Conversion Wash Buffer supplied with the 96-reaction kits (p/n 5271-0142) requires 32 mL of 100% ethanol.

CAUTION

For all washing steps in this section, do NOT spin the tubes in a centrifuge. Handle the tubes carefully to avoid any splashing.

- 1 Add 150 μ L of 1X Soft Conversion Wash Buffer to each tube of beads. Mix by pipetting up and down 15–20 times.

When handling multiple samples, Agilent recommends pouring 1X Soft Conversion Wash Buffer into a reagent reservoir and pipetting it using a multichannel pipette.

- 2 **Transfer the contents of each tube to a fresh strip tube**, then place the fresh strip tube on the magnet for 30 seconds or until the solution becomes clear (up to 1 minute). Remove and discard the supernatant from each tube, then remove the tubes from the magnet.
- 3 Add 150 μ L of 1X Soft Conversion Wash Buffer to each tube of beads. Mix by pipetting up and down 15–20 times.
- 4 Place the tubes on the magnet for 30 seconds or until the solution becomes clear (up to 1 minute). Remove and discard the supernatant from each tube, then use a 20 μ L pipette to remove any residual supernatant.
- 5 Remove the tubes from the magnet and proceed immediately to [“Step 4. DNA Repair.”](#)

Step 4. DNA Repair

- 1 Add 15 μL of freshly prepared 0.1X Soft Conversion Elution Buffer (prepared on [page 36](#)) to each tube. With caps closed, gently flick the tubes to ensure the beads go into solution (mixing by pipetting is not needed). Spin for 1 to 2 seconds to collect the liquid at the bottom of the tubes, making sure the beads do not settle.

NOTE

When handling multiple samples, Agilent recommends aliquoting 18 μL of 0.1X Soft Conversion Elution Buffer into the tubes of a fresh strip tube (one tube per sample). Then, use a multichannel pipette to transfer 15 μL of the buffer from the strip tube to the beads bound with converted library. This technique helps prevent the beads from drying out.

- 2 Program the thermal cycler to run the DNA elution thermal cycler program shown in [Table 25](#). Set the heated lid to 75°C. Load the strip tubes and run the program.

Table 25 Thermal cycler program for DNA elution

Step	Temperature	Time
Step 1	50°C	10 minutes
Step 2	10°C	Hold

- 3 Once the thermal cycler program for DNA elution reaches the 10°C hold step, remove the strip tubes from the thermal cycler. Add 5 μL of Soft Conversion Repair Solution to each tube (mixing is not needed). Spin for 1 to 2 seconds.
- 4 Place the tubes on the magnet for 30 seconds or until the solution becomes clear (up to 1 minute).
- 5 **Transfer all 20 μL of supernatant from each tube of the strip tube (carefully avoiding any bead carryover) to a fresh strip tube.** Proceed to [Chapter 6](#), "Indexing PCR, Library Purification, and Quality Assessment."

6

Indexing PCR, Library Purification, and Quality Assessment

Step 1. Indexing PCR [40](#)

Step 2. Library Purification [42](#)

Step 3. Library Quality Assessment and Quantification [43](#)

This chapter describes the steps for the indexing PCR, library purification, and quality assessment section of the Avida Methyl workflow. The captured libraries are amplified and indexed, and the indexed libraries are then purified with AMPure XP Beads. Finally, the purified, indexed libraries are quantified and analyzed for quality.

This chapter uses the reagents listed in [Table 26](#).

Table 26 Reagents for indexing PCR, library purification, and quality assessment

Reagent	Usage Notes	Kit/Storage Location
2X Methyl Amplification Mastermix (tube with black cap or bottle)	Thaw on ice.	Avida Methyl Reagent Box 1, stored at –20°C
Avida Index Primer Pairs	Primer pairs are provided in either a strip or a plate. Thaw strip or plate on ice.	Stored at –20°C
Nuclease-free Water (tube with clear cap or bottle)	—	Avida Methyl Reagent Box 2, stored at room temperature
AMPure XP Beads	Equilibrate at room temperature for 30 minutes.	Not provided
100% ethanol	—	Not provided
1X Low TE Buffer	—	Not provided
Qubit dsDNA HS or BR Assay Kit (optional)	—	Not provided
Nucleic acid analysis kit	—	Not provided

Step 1. Indexing PCR

CAUTION

Take precautions to avoid amplicon contamination during setup of the indexing PCR reactions. Review “[Important practices for preventing contamination](#)” on page 10.

- 1 Thaw the 2X Methyl Amplification Mastermix and Avida Index Primer Pairs on ice before use. Once thawed, gently vortex the 2X Methyl Amplification Mastermix at low speed, and briefly centrifuge the strip or plate of Avida Index Primer Pairs.
- 2 To each tube containing a converted library (from [step 5](#) on [page 38](#)), add 25 μL of 2X Methyl Amplification Mastermix to each tube. Mix by pipetting up and down 15–20 times. Briefly spin.
- 3 Add 5 μL of the appropriate Avida Index Primer Pair to each tube (use a pipette tip to pierce the seal of the tube or well containing the primer pair). Mix by pipetting up and down 15–20 times. Briefly spin.

Use a different index primer pair for each sample to be sequenced in the same lane.

The tubes now contain all reagents needed for indexing PCR. [Table 27](#) summarizes the contents in each PCR reaction.

Table 27 Indexing PCR reaction components

Reagent	Volume per PCR reaction
Converted library (from step 5 on page 38)	20 μL
2X Methyl Amplification Mastermix	25 μL
Avida Index Primer Pair	5 μL
Total	50 μL

NOTE

Index primer pairs are designed to be used in groups of 8. That is, pairs 1–8 are compatible with each other, as are pairs 9–16, 17–24, etc. For samples to be sequenced in the same lane, the best practice is to use primer pairs within a compatible set (e.g., 1–8). If you use primer pairs from across different compatibility sets, then you need to confirm that the selected pairs meet the compatibility requirements for Illumina sequencing. Specifically, make sure that none of the 16 nucleotide positions (8 on forward primer and 8 on reverse primer) contain only A and C nucleotides or only T and G nucleotides across all indexes in the set.

- 4 Program the thermal cycler to run the indexing PCR thermal cycler program shown in [Table 28](#). Set the heated lid to 103°C. Load the strip tubes and run the program.

Table 28 Thermal cycler program for indexing PCR

Step	Number of Cycles	Temperature	Time
Initial denaturation	1	98°C	45 seconds
Amplification stage 1	5	98°C	10 seconds
		62°C	30 seconds
		65°C	1 minute
Amplification stage 2	Variable, see Table 29 for recommendations	98°C	10 seconds
		65°C	1 minute
Final extension	1	65°C	1 minute
Final hold	1	4°C	Hold

Table 29 Indexing PCR cycle number recommendations for Amplification stage 2

Probes size of panel	Number of cycles based on 10 ng cfDNA input -- See Note below for other sample types/quantities
>100 kb	18 or fewer • Avida Methyl 3400 DMR Cancer Panel: 16 cycles
>50 kb to 100 kb	19
10 kb to 50 kb	20
<10 kb	21

NOTE

The cycle number recommendations in [Table 29](#) are based on a sample input of 10 ng of cfDNA. If using other sample types or input quantities, you may need to optimize the number of cycles using the points below for guidance.

- High-quality gDNA samples typically require 1–2 cycles more than cfDNA samples of the same input quantity.
- Low-quality gDNA samples (e.g., FFPE-derived gDNA) may require more cycles than high-quality gDNA samples. The number of additional cycles depends on the degree of sample degradation, with greater degradation necessitating more cycles.
- A sample input >50 ng typically requires 1–2 fewer cycles than a sample input of 10 ng.

- 5 Once the thermal cycler program for indexing PCR completes, remove the strip tubes containing the amplified, indexed libraries from the thermal cycler. Put on ice.

Stopping Point If you do not continue to the next step, store the indexed libraries at 4°C overnight or at –20°C for up to 72 hours. Avoid prolonged storage >72 hours.

Step 2. Library Purification

- 1 Equilibrate the AMPure XP Beads at room temperature for 30 minutes before use.
- 2 Prepare 400 μL of fresh 80% ethanol (dilute 100% ethanol with Nuclease-free Water) per library, multiplied by the number of libraries plus a 10% overage.
- 3 Thoroughly vortex the stock of AMPure XP Beads to mix. Add 50 μL (1X volume) of the AMPure XP Beads to each tube containing an amplified, indexed library. Mix by pipetting up and down 15–20 times.
- 4 Incubate for 5 minutes at room temperature.
- 5 Place the strip tubes on a magnet for 2 minutes or until the solution is clear. Remove and discard the supernatant from each tube.
- 6 With the tubes still on the magnet, add 200 μL of freshly prepared 80% ethanol to each tube without disturbing the beads. Incubate for 30 seconds, then carefully remove and discard the supernatant from each tube.
- 7 Perform an additional 80% ethanol wash by repeating [step 6](#) for a total of 2 washes.
- 8 Use a 20 μL pipette to remove all residual 80% ethanol. Air dry the beads for up to 3 minutes, keeping the tubes on the magnet with the lids off. Do not over-dry the beads.
- 9 Remove the strip tubes from the magnet, then add 23 μL of 1X Low TE Buffer to each tube. Mix by pipetting up and down 15–20 times. Briefly spin.
- 10 Incubate at room temperature for 2 minutes.
- 11 Place the strip tubes on a magnet for 2 minutes or until the solution becomes clear.
- 12 **Transfer 20 μL of the eluate from each tube to a new tube**, being careful to avoid bead carryover.

Stopping Point If you do not continue to the next step, store the indexed libraries at 4°C overnight or at –20°C for prolonged storage.

Step 3. Library Quality Assessment and Quantification

- 1 (Optional) Measure the libraries using a fluorescence-based method for DNA quantification, such as the Qubit dsDNA HS Assay Kit or Qubit dsDNA BR Assay Kit. Follow the manufacturer's instructions for the instrument and assay kit.
- 2 Analyze a sample of each library using one of the platforms listed in [Table 30](#). Follow the instructions in the linked user guide provided for each assay.

Table 30 Library analysis options

Analysis platform	Assay used at this step (links go to assay instructions)	Amount of library sample to analyze
Agilent 4200/4150 TapeStation system	D1000 ScreenTape	1 µL of sample mixed with 3 µL of D1000 sample buffer*
Agilent 5200, 5300, or 5400 Fragment Analyzer system	One of the following: NGS Fragment Kit (1-6000 bp) Small Fragment Kit (1-1500 bp) HS NGS Fragment Kit (1-6000 bp) HS Small Fragment Kit (1-1500 bp)	2 µL of sample

* If the sample input for the workflow was near the top of the input range of 100 ng, you may need to further dilute the sample prior to analysis to make sure the quantity is in the dynamic range for the TapeStation system.

Each analysis method provides an electropherogram showing the size distribution of fragments in the sample and tools for determining the concentration of DNA in the sample. See [Table 31](#) for fragment size distribution guidelines.

To illustrate typical results, representative electropherograms generated using the TapeStation system are provided in [Figure 2](#) (cfDNA sample) and [Figure 3](#) (gDNA sample). Representative electropherograms generated on the Fragment Analyzer system are provided in [Figure 4](#) (cfDNA sample) and [Figure 5](#) (gDNA sample).

Determine the concentration of the library DNA by integrating under the entire peak.

Table 31 Library qualification guidelines

Input DNA type	Expected DNA fragment size peak position
cfDNA	320 bp (see Figure 2 and Figure 4 for sample electropherograms)
Fragmented gDNA*	300–350 bp (see Figure 3 and Figure 5 for sample electropherograms)

* Refers to high-quality gDNA samples. For gDNA derived from FFPE samples, the fragment sizes may be smaller.

Stopping Point

If you do not sequence the libraries immediately, store the tubes at 4°C overnight or at –20°C for prolonged storage.

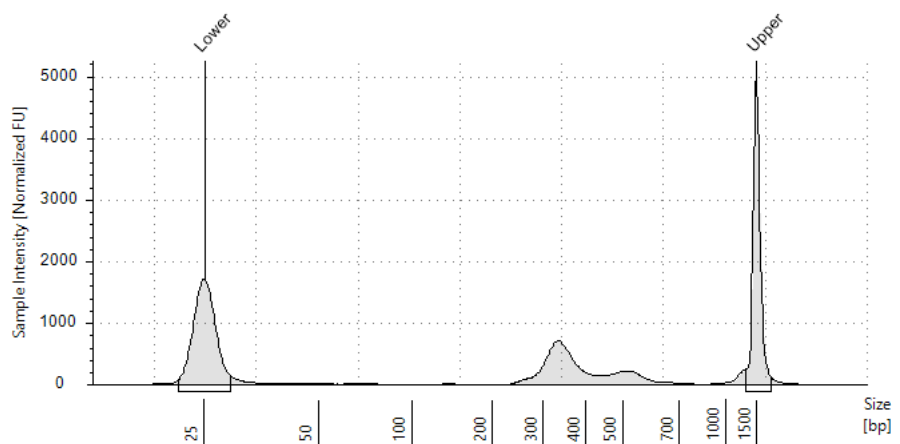


Figure 2 Library prepared from 10 ng of a cfDNA sample hybridized with the Avida Methyl 3400 DMR Cancer Panel and analyzed using a D1000 ScreenTape assay on the TapeStation.

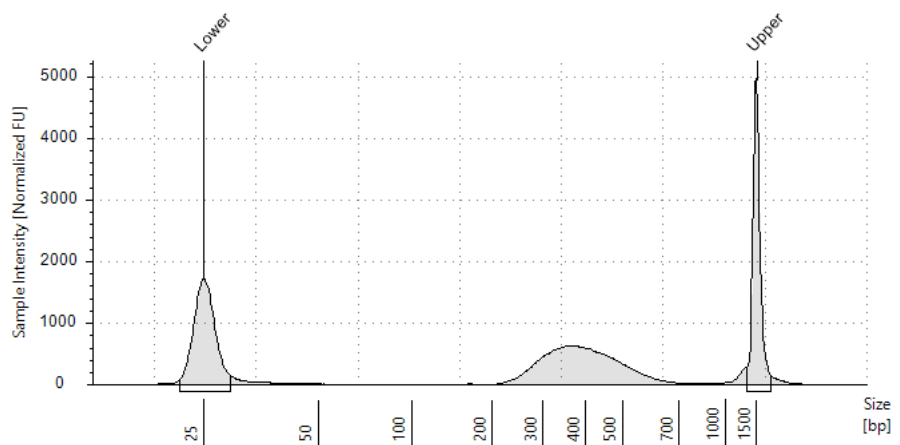


Figure 3 Library prepared from 10 ng of a high-quality gDNA sample hybridized with the Avida Methyl 3400 DMR Cancer Panel and analyzed using a D1000 ScreenTape assay on the TapeStation.

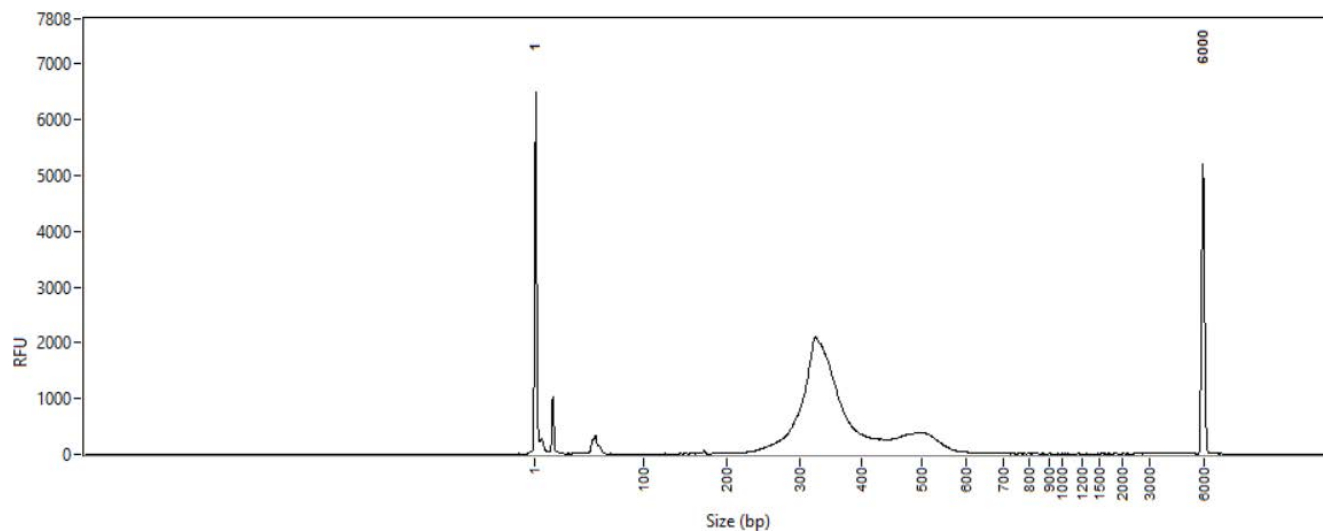


Figure 4 Library prepared from 10 ng of a cfDNA sample hybridized with the Avida Methyl 3400 DMR Cancer Panel and analyzed using the NGS Fragment Kit on the Fragment Analyzer.

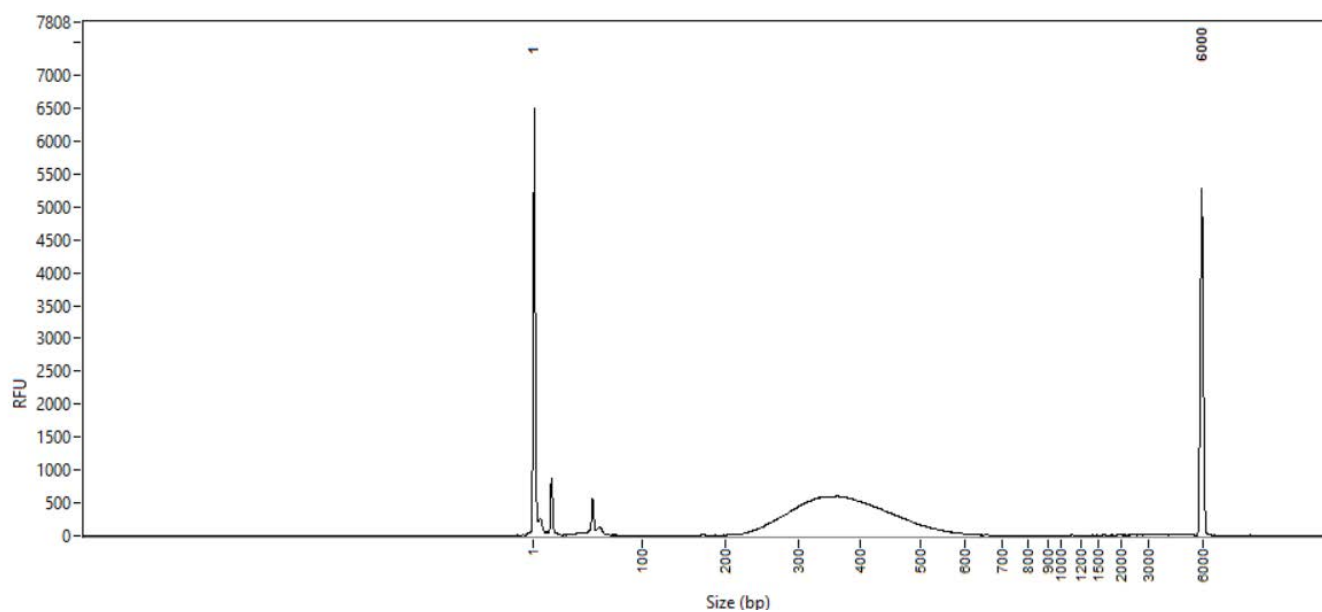


Figure 5 Library prepared from 10 ng of a high-quality gDNA sample hybridized with the Avida Methyl 3400 DMR Cancer Panel and analyzed using the NGS Fragment Kit on the Fragment Analyzer.

7 Sequencing and NGS Analysis

Step 1. Pool samples for multiplexed sequencing [47](#)

Step 2. Prepare the sequencing samples [48](#)

Step 3. Sequence the libraries [49](#)

Step 4. Process and analyze the reads [51](#)

This chapter contains guidance on library sequencing and analysis. Refer to your specific Illumina sequencer's user guide for specific instructions on how to perform sequencing.

Step 1. Pool samples for multiplexed sequencing

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

Combine the libraries such that each indexed library is present in equimolar amounts in the pool using one of the following methods. Use the diluent specified by your sequencing provider, such as 1X Low TE Buffer, for the dilution steps.

Method 1: Dilute each library 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 pool.

Method 2: Starting with library 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 1X Low TE Buffer. The formula below is provided for determination of the amount of each indexed sample to add to the pool.

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

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

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

$\#$ is the number of indexes, and

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

Table 32 shows an example of the amount of 4 index-tagged samples (of different concentrations) and 1X Low TE Buffer needed for a final volume of 20 µL at 10 nM DNA.

Table 32 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
1X Low TE Buffer					7.6

Step 2. Prepare the sequencing samples

The final Avida library pool is ready for sequencing using standard Illumina paired-end primers and chemistry. Each fragment in the prepared library contains one target insert surrounded by sequence motifs required for multiplexed sequencing using the Illumina platform ([Figure 6](#)).

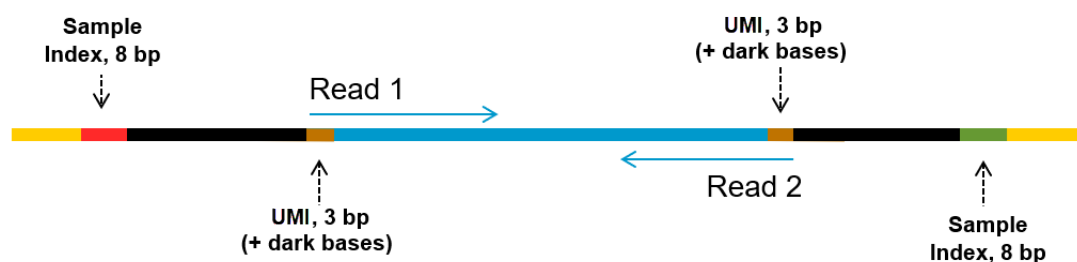


Figure 6 Content of Avida sequencing library. Each fragment contains one target insert (blue) surrounded by the Illumina paired-end sequencing elements (black), unique dual sample indexes (red and green), duplex UMIs (brown), and the library PCR primers (yellow).

Proceed to cluster amplification using the appropriate Illumina Paired-End Cluster Generation Kit and sequence the libraries using an Illumina instrument. Consult Illumina's documentation for kit configuration and seeding concentration guidelines.

Seeding concentration and cluster density may need to be optimized based on the DNA fragment size range for the library and on the desired output and data quality. Begin optimization using a seeding concentration in the middle of the range provided by Illumina. Follow Illumina's recommendation for a PhiX control in a low-concentration spike-in for improved sequencing quality control.

Step 3. Sequence the libraries

Proceed to cluster amplification using the appropriate Illumina Paired-End Cluster Generation Kit. [Table 33](#) provides guidelines for use of several Illumina instrument and chemistry combinations suitable for processing Avida target-enriched NGS libraries. For other Illumina NGS platforms, consult Illumina's documentation for kit configuration and seeding concentration guidelines.

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

Table 33 Illumina Kit Configuration Selection Guidelines

Instrument	Run Type	Read Length	SBS Kit Configuration	Chemistry	Seeding Concentration
MiSeq	All Runs	2 × 100 bp or 2 × 150 bp	300 Cycle Kit	v2	9–10 pM
			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

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 34](#) showing example settings for 2 × 150 bp sequencing.

Table 34 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. See [“Online resource for Avida index sequences”](#) on page 13 for information on downloading the Avida index sequences spreadsheet.
- No custom primers are used for Avida library sequencing. Leave all Custom Primers options for Read 1, Read 2, Index 1 and Index 2 primers cleared/deselected during run setup.
- 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 adaptors, including the UMIs in the adaptor sequences.
- 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 Avida index sequences provided online 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 Avida run setup in your selected application (e.g., generating index files or Sample Sheet templates), contact the Agilent Technical Support team (see [page 2](#)) or your local representative.

Step 4. Process and analyze the reads

Typical sequencing data processing steps are outlined below.

- 1 Generate demultiplexed FASTQ files for each sample using Illumina's bcl2fastq, BCL Convert, or a similar software application. This process generates paired-end reads based on the dual indexes and removes sequences with incorrectly paired P5 and P7 indexes. Do not use the UMI trimming options offered in Illumina's demultiplexing software.

NOTE

If your sequence analysis pipeline excludes UMIs, you can remove the first 5 bases from Read 1 and Read 2 by masking or trimming before proceeding to downstream analysis.

If demultiplexing using bcl2fastq, UMIs may be masked by including the base mask **N5Y*,I8,I8,N5Y*** (where * is replaced with the remaining read length after subtracting the 5 masked bases, e.g., use **N5Y146,I8,I8,N5Y146** for 2x150 NGS set up as shown in [Table 34](#) on page 49). The sum of the values following N and Y must match the read length value in the RunInfo.xml file.

If demultiplexing using BCL Convert, UMIs may be trimmed by including the following string in the sample sheet header: **OverrideCycles,N5Y*,I8,I8,N5Y*** (where * is replaced with read length after trimming, e.g., use **N5Y146,I8,I8,N5Y146** for 2x150 NGS set up as shown in [Table 34](#) on page 49). The sum of the values following N and Y must match the read length value in the RunInfo.xml file.

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

- 2 Use a suitable processing tool of your choice to trim and collect inline UMIs from each sequencing read.
 - Inline UMIs are added to both ends of the DNA inserts in the assay. To collect the UMIs and process them for analysis, one approach is to trim off 5 bases at the beginning of each read, take the first 3 of the 5 bases as UMI, and discard the remaining 2.
- 3 (Optional) Perform additional adaptor trimming with a tool such as Trimmomatic or Cutadapt, or hard trim both the R1 and R2 reads on the 3' end to approximately 100PE. This additional trimming may help improve the alignment rate, especially for DNA inserts that are shorter than 200 bp.
- 4 Use Bismark or other suitable tool to align the paired end reads and generate a BAM file. An example command is below.

```
[path_to_bismark_command] --bowtie2 \
  [path_to_bismark_converted_reference_genome] \
  [path_to_samtools] \
  -1 R1.fq -2 R2.fq \
  --nucleotide_coverage \
  --output_dir [path_to_output_dir] \
  --temp_dir [path_to_output_dir] --basename bs_aln
```

- 5 Remove duplicate reads based on strand, genomic location, and, optionally, UMI. For UMI based deduplication, use a tool such as Umi-Grinder. An example command is below.

```
[path_to_Umi-Grinder]/UmiBam -p --bam --umi --mm 1 \  
--samtools_path [path_to_samtools] [aligned_reads].bam
```
- 6 (Optional) Analyze deduplicated reads to form duplex UMI consensus sequences.
- 7 Collect QC metrics, such as raw coverage, deduplicated coverage, alignment percentage, and duplex rate.
- 8 Extract CpG (5'C-phosphate-G-3') and CHX methylation information from the deduplicated reads with a tool such as *bismark_methylation_extractor*.

For CHX, H is A, C or T, while X can be A, C, G or T.
- 9 (Optional) Perform methylation profiling of CpGs in the panel regions.
- 10 (Optional) Generate a report containing the QC metrics and methylation profile.

8 Reference

Reagent Kit Contents	54
Avida Index Primer Pair Information	56
Avida Custom AddOn Panels	58
Troubleshooting Guide	59
Quick Reference Protocol	60

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

Reagent Kit Contents

[Table 2](#) on page 12 lists the Avida Methyl reagent kits and their sub-kit boxes. Detailed contents of those sub-kit boxes are shown in [Table 35](#) through [Table 38](#).

Avida Sub-Kit Details

Table 35 Avida Methyl Reagent Box 1

Kit Component	16 Reaction Kit (p/n 5282-0145)	96 Reaction Kit (p/n 5282-0146)
End Prep Buffer	tube with purple cap	tube with purple cap
End Prep Enzyme	tube with blue cap	tube with blue cap
Ligation Buffer	tube with green cap	bottle
Ligation Enzyme	tube with yellow cap	tube with yellow cap
Adaptor for ILM	tube with orange cap	tube with orange cap
Hyb Blocker	tube with red cap	tube with red cap
Hyb Buffer	tube with clear cap	bottle
Hyb Enhancer	amber tube with green cap	tube with green cap
Soft Conversion Reagent A	amber tube with amber cap	tube with amber cap
2X Methyl Amplification Mastermix	tube with black cap	bottle

Table 36 Avida Methyl Reagent Box 2

Kit Component	16 Reaction Kit (p/n 5282-0147)	96 Reaction Kit (p/n 5282-0148)
Library Wash Buffer	bottle	bottle
Hyb Wash Buffer 1	bottle	bottle
Hyb Wash Buffer 2	bottle	bottle
Nuclease-Free Water	tube with clear cap	bottle

Table 37 Avida Methyl and Duo Reagent Box 3

Kit Component	16 Reaction Kit (p/n 5282-0149)	96 Reaction Kit (p/n 5282-0150)
Soft Conversion Binding Beads	bottle	bottle
5X Soft Conversion Wash Buffer	bottle	bottle
Soft Conversion Elution Buffer	tube with green cap	tube with green cap
Soft Conversion Repair Solution	tube with yellow cap	tube with yellow cap
Nuclease-Free Water	tube with clear cap	bottle

Reference

Avida Sub-Kit Details

Table 38 Soft Conversion Reagent B

Kit Component	16 Reaction Kit (p/n 5280-0188)	96 Reaction Kit (p/n 5280-0189)
Soft Conversion Reagent B*	amber tube with blue cap	tube with blue cap

* Soft Conversion Reagent B may be shipped as part of the Avida Methyl and Duo Reagent Box 3 or it may be shipped in its own separate bag at room temperature.

Table 39 Avida Beads Box

Kit Component	16 Reaction Kit (p/n 5282-0143)	96 Reaction Kit (p/n 5282-0144)
Library Binding Beads	tube with white cap	bottle
Capture Beads	tube with amber cap	tube with amber cap

Avida Index Primer Pair Information

The Avida Index Primer Pairs are provided pre-combined in the well of either a strip tube or a plate (see [Table 40](#)). Each member of the primer pair contains a unique 8-bp P5 or P7 index, resulting in dual-indexed NGS libraries. One primer pair (forward and reverse primers) is provided as a single-use aliquot in each well of an 8-well strip tubes or 96-well plate.

See [“Online resource for Avida index sequences”](#) on page 13 for information on downloading the Avida index sequences spreadsheet.

Table 40 Avida Index Primer Pairs for ILM Content

Kit Component	16 Reaction Kit Format	96 Reaction Kit Format
Avida Index Primer Pairs for ILM	Black 8-well strip tube with index pairs 1-8, AND Blue 8-well strip tube with index pairs 9-16 -- OR -- Red 8-well strip tube with index pairs 17-24, AND White 8-well strip tube with index pairs 25-32	Clear 96-well plate with index pairs 1–96 -- OR -- Blue 96-well plate with index pairs 97–192

Index Primer Pair Strip Tube and Plate Maps

- The black strip contains Index Primer Pairs 1-8, with pair #1 supplied in the well proximal to the numeral **1** etched on the strip’s plastic end tab.
- The blue strip contains Index Primer Pairs 9-16, with pair #9 supplied in the well proximal to the numeral **9** etched on the strip’s plastic end tab.
- The red strip contains Index Primer Pairs 17-24, with pair #17 supplied in the well proximal to the numeral **17** etched on the strip’s plastic end tab.
- The white strip contains Index Primer Pairs 25-32, with pair #25 supplied in the well proximal to the numeral **25** etched on the strip’s plastic end tab.

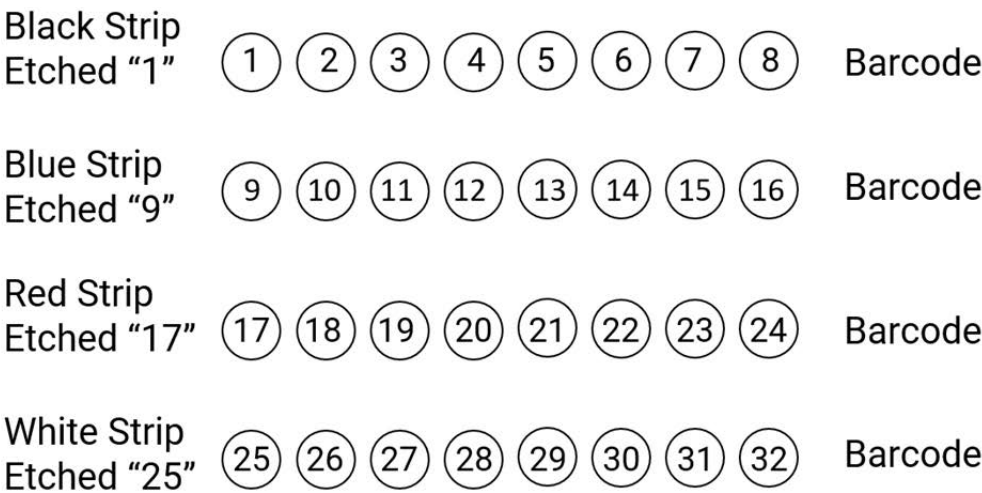


Figure 7 Map of the Avida Index Primer Pairs strip tubes provided with 16 reaction kits

[Table 41](#) and [Table 42](#) show the plate positions of the Avida Index Primer Pairs provided with 96 reaction kits.

Table 41 Plate map for Avida Index Primer Pairs 1-96, provided in clear plate

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

Table 42 Plate map for Avida Index Primer Pairs 97-192, provided in blue plate

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

Avida Custom AddOn Panels

When including an Avida Methyl Custom AddOn Panel in the hybridization, make the protocol adjustments described below.

CAUTION

Make sure that the Custom AddOn panel is compatible with the Avida catalog panel you are using in the workflow. If needed, check with Agilent [Technical Support](#) or your local sales representative for design ID compatibility. Custom AddOn panels are limited to 25 kb in size.

During preparation of the Hyb Mix 1 on [page 25](#), substitute the reagents and volumes in [Table 15](#) with those shown in [Table 43](#) below.

Table 43 Preparation of Hyb Mix 1 using Avida Methyl Custom AddOn Panel

Reagent	Volume for 1 reaction	Volume for 8 reactions (includes excess)	Volume for 16 reactions (includes excess)
Nuclease-Free Water	13.5 µL	121.5 µL	243 µL
Avida Methyl 3400 DMR Cancer Panel (catalog)	4 µL	36 µL	72 µL
Avida Methyl Custom AddOn Panel	4 µL	36 µL	72 µL
Hyb Blocker	2.5 µL	22.5 µL	45 µL
Total	24 µL	216 µL	432 µL

Troubleshooting Guide

If yield of libraries is lower than previous captures

- Make sure that the Library Binding Beads (used in [“Step 3. Library bead binding”](#) on page 25) have equilibrated to room temperature before use.
- During [“Step 1. Indexing PCR”](#) on page 40, the number of cycles used in for Amplification stage 2 may require optimization. Refer to the Note on [page 41](#) for guidelines.
- Take care throughout the workflow to minimize sample loss during pipetting and transfer steps.
- When mixing samples with beads, make sure the reactions are well mixed.

If peak positions in the electropherogram are not as expected

- The indexing PCR reactions may have been contaminated with amplicon from another assay. Maintain separate work areas for pre-PCR and post-PCR steps. If maintaining separate areas is not an option, clean the at-risk surfaces with 10% bleach, then wipe down with water to remove bleach residue. Review [“Important practices for preventing contamination”](#) on page 10.

If low percent on-target is observed in library sequencing results

- During [“Step 3. Library bead binding”](#) on page 25, include the optional second wash (see [step 10](#) on [page 26](#)) and make sure to remove all residual Library Wash Buffer during the washes. The following techniques can be used to collect residual supernatant at the bottom of the tubes:
For magnets that collect beads on tube sides: Tap the magnet stand on the bench 5 times.
For magnets that collect beads at tube bottoms: Remove tubes from magnet and briefly spin. Return tubes to magnet before removing liquid.
- Following hybridization, make sure that the Capture Beads are fully resuspended during the wash steps with Hyb Wash Buffer 1 and Hyb Wash Buffer 2 and make sure to remove all residual wash buffer (see [“Step 4. Washing \(Bead-Bound DNA Hybrids\)”](#) on page 32). Refer to techniques above for collecting residual supernatant based on magnet type.

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

- Make sure that the hybridization temperature used in the hybridization thermal cycle program is 60°C as directed in [Table 18](#) on page 28. Additionally, make sure that the thermal cycler performs the washing incubation at 60°C as directed in [Table 20](#) on page 31.

If conversion rate is low

- During [“Step 1. Soft Conversion”](#) on page 34, all residual supernatant must be removed from the beads before addition of the Soft Conversion Elution Buffer. The following techniques can be used to collect residual supernatant at the bottom of the tubes:
For magnets that collect beads on tube sides: Tap the magnet stand on the bench 5 times.
For magnets that collect beads at tube bottoms: Remove tubes from magnet and briefly spin. Return tubes to magnet before removing liquid.
- During [“Step 1. Soft Conversion”](#) on page 34, the samples need to be well mixed with Soft Conversion Mix. Make sure to pipette the mixture up and down at least 15–20 times.
- The thermal cycler program for the soft conversion reactions is provided in [Table 23](#) on page 35. Use the durations and temperatures as listed in the table.
- During [“Step 3. Washing \(Bead-Bound Converted Libraries\)”](#) on page 37, make sure that all residual 1X Soft Conversion Wash Buffer is removed from the bead pellet before adding the 0.1X Soft Conversion Elution Buffer. Refer to techniques above for collecting residual supernatant based on magnet type.

Quick Reference Protocol

An abbreviated summary of the protocol steps is provided below for experienced users. Use the complete protocol until you are familiar with all of the protocol details such as reagent mixing instructions and instrument settings.

A new print-ready, graphics-based quick start protocol is available on the Agilent website:
<https://www.agilent.com/cs/library/usermanuals/public/G9419-90500.pdf>.

Step	Summary of Conditions
<i>Before starting a protocol:</i> Pre-heat Hyb Wash Buffer 1	Heat stock bottle to 50°C in water bath or heat block In a 50°C water bath: Heat for 10 minutes, then transfer the bottle to a 50°C heat block until use On a 50°C heat block: Leave the bottle on the 50° heat block until use
PCR-free Library Prep	
Prepare End Prep Mix	Per 8 reactions: 63 µL End Prep Buffer + 27 µL End Prep Enzyme Per 16 reactions: 126 µL End Prep Buffer + 54 µL End Prep Enzyme Keep on ice or use immediately
End-Repair and dA-Tail the DNA fragments	50 µL DNA (cfDNA or fragmented gDNA) + 10 µL End Prep Mix Mix and spin, then keep on ice Incubate in thermal cycler (heated lid 75°C): 30 min @ 20°C, 30 min @ 65°C, Hold @ 4°C
Prepare Ligation master mix	Per 8 reactions: 225 µL Ligation Buffer + 54 µL Ligation Enzyme Per 16 reactions: 450 µL Ligation Buffer + 108 µL Ligation Enzyme
Ligate adaptor	Add 5 µL Adaptor for ILM to each sample Then, add 31 µL Ligation Master Mix Mix and spin Incubate in thermal cycler (heated lid off): 30 min @ 20°C, Hold @ 4°C
Prepare reagents	Library Binding Beads: Room temperature for at least 15 minutes Hyb Blocker and Hyb Enhancer: Thaw then keep at room temperature Avida Methyl Panel: Thaw then keep on ice Hyb Buffer: 37°C for at least 10–20 minutes, then keep at room temperature
Bind to Library Binding Beads	Vortex Library Binding Beads Add 87 µL Library Binding Beads to each sample Mix and spin without letting beads settle Incubate 10 min @ room temperature
<i>During 10-min incubation:</i> Prepare Hyb Mix 1	Per 8 reactions: 157.5 µL Nuclease-free Water + 36 µL Avida Panel + 22.5 µL Hyb Blocker Per 16 reactions: 315 µL Nuclease-free Water + 72 µL Avida Panel + 45 µL Hyb Blocker Keep at room temperature
Wash	After 10-min incubation, place tubes on magnet, discard supernatant Wash beads in 180 µL Library Wash Buffer Place tubes on magnet, discard supernatant (Optional) Repeat wash with Library Wash Buffer Remove ALL remaining supernatant Remove samples from magnet
Target Capture	
Add Hyb Mix 1	Add 24 µL Hyb Mix 1 Mix and spin

Step	Summary of Conditions
Prepare Hyb Mix 2	Per 8 reactions: 270 μ L Hyb Buffer + 54 μ L Hyb Enhancer Per 16 reactions: 540 μ L Hyb Buffer + 108 μ L Hyb Enhancer
Hybridize	Add 36 μ L Hyb Mix 2 Mix and spin without letting beads settle Incubate in thermal cycler (heated lid 103°C): 2 min @ 98°C, $\geq 2.5^\circ\text{C}/\text{second}$ ramp down to 60°C, 60 min @ 60°C, Hold @ 4°C
<i>During hybridization program:</i> Prepare Capture Beads	Mix Capture Beads well Calculate volume of Capture Beads needed (8 μ L/sample + overage) Transfer that volume into 1.5-mL tube Place tube on magnet, discard supernatant, then remove tube from magnet Wash beads 2 \times in 100 μ L preheated Hyb Wash Buffer 1, discard supernatant after each wash (If processing >16 reactions, scale the volume of Hyb Wash Buffer 1 to use 100 μ L per 16 reactions; see Table 19 on page 29) Remove tube from magnet Resuspend Capture Beads in Hyb Buffer at original volume calculated above, and mix Add 8 μ L resuspended Capture Beads to each tube of a fresh strip tube
Bead capture	After hybridization, spin tubes and place on magnet Transfer each supernatant to tubes containing 8 μ L-aliquots of prepared Capture Beads Mix and spin without letting beads settle Incubate in thermal cycler held @60°C (heated lid 75°C) After 10 minutes @ 60°C, remove tubes but leave thermal cycler running Place tubes on magnet, discard supernatant Remove samples from magnet
Wash	Wash beads in 150 μ L preheated Hyb Wash Buffer 1 Place tubes on magnet and discard supernatant Remove samples from magnet Add 100 μ L Hyb Wash Buffer 1 and mix Transfer each sample to a fresh strip tube and cap each tube Incubate in thermal cycler held @60°C (heated lid 75°C) After 3 minutes, place on magnet, discard supernatant Remove samples from magnet Add 150 μ L Hyb Wash Buffer 2, and mix
Soft Conversion and Repair	
Prep samples for conversion	Briefly spin tubes Place tubes on magnet and discard supernatant Remove ALL remaining supernatant Remove samples from magnet Add 11.5 μ L Soft Conversion Elution Buffer, and pulse vortex to mix Incubate 10 min @ room temperature
<i>During 10-min incubation:</i> Prepare Soft Conversion Mix	Make sure Soft Conversion Reagent A and Soft Conversion Reagent B are well mixed Per 8 reactions: 31.5 μ L Soft Conversion Reagent A + 180 μ L Soft Conversion Reagent B Per 16 reactions: 63 μ L Soft Conversion Reagent A + 360 μ L Soft Conversion Reagent B Mix and spin
Perform soft conversion	After the 10-minute incubation, place tubes on magnet Transfer supernatant to fresh tubes Add 23.5 μ L of freshly prepared Soft Conversion Mix Mix and spin Incubate in thermal cycler (heated lid 85°C): 25 min @ 75°C, Hold @ 10°C Proceed to next step as soon as thermal cycler reaches 10°C hold

Step	Summary of Conditions
Set up bead purification	Add 150 μ L of Soft Conversion Binding Beads Mix and spin without letting beads settle Incubate 10 min @ room temperature
<i>During 10-min incubation:</i> Prepare 0.1X Soft Conversion Elution Buffer	Per 8 reactions: 16.5 μ L Soft Conversion Elution Buffer + 148.5 μ L Nuclease-Free Water Per 16 reactions: 33 μ L Soft Conversion Elution Buffer + 297 μ L Nuclease-Free Water Mix and spin
Wash	After the 10-minute incubation, place tubes on magnet and discard supernatant Add 150 μ L of 1X Soft Conversion Wash Buffer and mix Transfer contents to fresh tubes Place tubes on magnet and discard supernatant Add 150 μ L of 1X Soft Conversion Wash Buffer and mix Place tubes on magnet and discard supernatant Remove tubes from magnet
Repair	Add 15 μ L of freshly prepared 0.1X Soft Conversion Elution Buffer Mix by flicking the tubes then spin without letting beads settle Incubate in thermal cycler (heated lid 75°C): 10 min @ 50°C, Hold @ 10°C As soon as thermal cycler reaches 10°C hold, add 5 μ L of Soft Conversion Repair Solution Spin (without mixing) Place tubes on magnet Transfer supernatant to fresh tubes
Indexing PCR, Library Purification, and Quality Assessment	
Set up PCR reactions	Add 25 μ L 2X Methyl Amplification Mastermix Add 5 μ L primer pair Mix and spin
Perform indexing PCR	Incubate in thermal cycler (heated lid 103°C): 45 sec @98°C 5 cycles: 10 sec @98°C, 30 sec @ 62°C, 1 minute @65°C Additional cycles (as described in Table 29): 10 sec @98°C, 1 min @65°C 1 min @65°C Hold @4°C
Purify libraries	Add 50 μ L of AMPure XP Beads and mix Incubate 5 min @ room temperature Place on magnet, discard supernatant, keep samples on magnet Wash beads 2x in 200 μ L 80% ethanol, discard supernatant after each wash Remove all residual ethanol then air dry the beads up to 3 minutes (on magnet, lids open) Remove samples from magnet Add 23 μ L of 1X Low TE Buffer Mix and spin Incubate at room temperature for 2 min then put samples back on magnet Transfer the eluate to fresh tube
Assess quality and quantity	(Optional) Quantification using fluorescence-based method (e.g., Qubit) Quality assessment on TapeStation or Fragment Analyzer Libraries are ready for sequencing or storage (4°C overnight or at -20°C for prolonged storage)

In This Book

This guide provides instructions for using Avida Methyl Reagent Kits for library preparation, target capture, and soft conversion.

