

 **BD Rhapsody™ System**
mRNA Whole Transcriptome Analysis
(WTA) and Sample Tag
Library Preparation Protocol

Copyrights

No part of this publication may be reproduced, transmitted, transcribed, stored in retrieval systems, or translated into any language or computer language, in any form or by any means: electronic, mechanical, magnetic, optical, chemical, manual, or otherwise, without prior written permission from BD.

The information in this guide is subject to change without notice. BD reserves the right to change its products and services at any time. Although this guide has been prepared with every precaution to ensure accuracy, BD assumes no liability for any errors or omissions, nor for any damages resulting from the application or use of this information. BD welcomes customer input on corrections and suggestions for improvement.

Patents and Trademarks

For US patents that may apply, see bd.com/patents.

BD, the BD Logo and BD Rhapsody are trademarks of Becton, Dickinson and Company or its affiliates. All other trademarks are the property of their respective owners. © 2024 BD. All rights reserved.

Regulatory information

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

History

Revision	Date	Change made
23-24119(01)	2021-12	Initial release.
23-24119(02)	2022-11	Updated for BD Rhapsody™ Enhanced Cell Capture Beads version 2.0. In the Appendix, added BD® Flex SMK sequences.
23-24119(03)	2024-11	Added the BD® OMICS-One Dual Index Kit and BD Rhapsody™ Enhanced Cartridge Reagent Kit v3. Added Workflows chapter. Updated Time Considerations workflow. Updated sequencing section. Added sequencing recommendation. Added BD Rhapsody™ oligo sequences to appendix.

Contents

Introduction	5
Symbols	5
Protocol kits	6
Workflows	7
WTA library amplification workflow	7
Sample Tag library amplification workflow	7
Required and recommended materials	8
Required reagents	8
Recommended consumables	8
Equipment	9
Best practices	9
Additional documentation	10
Safety information	10
Time considerations	11
Procedure	12
1. WTA library amplification	12
Before you begin	12
1.1 Random priming and extension (RPE)	13
1.2 RPE cleanup	16
1.3 Additional RPE cleanup for cell input <5,000 PBMC cells	17
1.4 RPE PCR	18
1.5 RPE PCR cleanup and quality check	19
1.6 WTA index PCR	22
1.7 WTA index PCR cleanup and quality check	23
1.8 Additional WTA index PCR cleanup	27
2. Sample Tag library amplification	28
2.1 Sample Tag PCR1	28
2.2 Sample Tag PCR1 cleanup	30
2.3 Sample Tag PCR2	31
2.4 Sample Tag PCR2 cleanup and quantify	31
2.5 Sample Tag index PCR	33
2.6 Sample Tag index PCR cleanup and quality check	35
Sequencing	37
Read requirements for libraries	37
Sequencing recommendations	38
Sequencing analysis pipeline	38
Troubleshooting	39
Library preparation	39
Sequencing	41
Appendix	42
Oligonucleotides in BD Rhapsody™ Whole Transcriptome Analysis Amplification Kit	42

Sample Tag sequences	44
BD® Flex SMK sequences	46
Contact Information	50

Introduction

This protocol provides instructions on creating single-cell whole transcriptome mRNA and Sample Tag libraries after cell capture on the BD Rhapsody™ HT Single-Cell Analysis System or the BD Rhapsody™ HT Xpress System for sequencing on various sequencers. For complete instrument procedures and safety information, refer to the *BD Rhapsody™ HT Single-Cell Analysis System Instrument User Guide* or the *BD Rhapsody™ HT Xpress System Instrument User Guide for Scanner-Free Workflow*.

The cDNA of mRNA and Sample Tag targets is first encoded on BD Rhapsody™ Enhanced Cell Capture Beads, as described in the instrument user guides. At the same time, the barcode information from BD Rhapsody™ Enhanced Cell Capture Beads is also added to Ab-Oligos during reverse transcription, which enables amplification of Sample Tags in solution. To generate the mRNA and Sample Tag sequencing libraries, the extended Sample Tags are first denatured from the BD Rhapsody™ Enhanced Cell Capture Beads, which are later amplified separately through a series of PCR steps. Meanwhile, the whole transcriptome amplification library is generated directly from the BD Rhapsody™ Enhanced Cell Capture Beads using a random priming approach, followed by an index polymerase chain reaction (PCR) step. The whole transcriptome mRNA and Sample Tag libraries can be combined together for sequencing on various sequencers.

This protocol is intended to provide a method to screen RNA expression of single cells using a 3' whole transcriptome analysis (WTA) approach through the BD Rhapsody™ WTA Amplification Kit for samples that have been labeled using Sample Tags from the BD® Single Cell Multiplexing Kit. The data set generated from this protocol can be used to generate a custom panel for subsequent 3' Targeted mRNA sequencing. Specifically, the protocol outlines how to generate whole transcriptome libraries for cell-capture inputs between 100 to 100,000 resting peripheral blood mononuclear cells (PBMCs) per sample for library generation. For cell-capture inputs between 100 to <5,000 cells per sample, there are sections in the protocol for additional cleanups. For cell types other than resting PBMCs, protocol optimization might be required by the user.

Symbols

The following symbols are used in this guide:

Symbol	Description
	Important information for maintaining measurement accuracy or data integrity.
	Noteworthy information.
	Procedural stopping point.

Protocol kits

Before you begin, ensure that you have the correct kits for this protocol. Matching cap colors indicate you have the correct kit, along with the catalog numbers found in the [Required and recommended materials \(page 8\)](#) section.

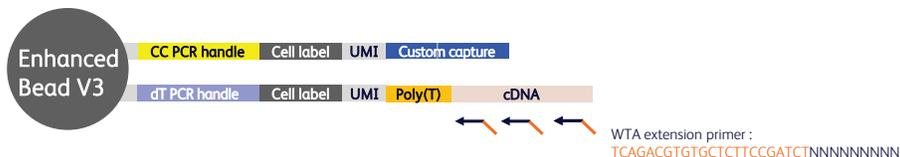
BD Rhapsody™ WTA Amplification Kit (-20 °C)		
Cap Color	Name	Quantity
●	WTA extension primers	1
●	WTA extension buffer	1
●	WTA extension enzyme	1
●	10 mM dNTP	1
○	Nuclease-free water	3
●	Bead RT/PCR enhancer	1
○	WTA amplification primer	1
○	PCR master mix	1
○	Universal oligo	2
●	Sample Tag PCR1 primer	2
●	Sample Tag PCR2 primer	1
●	BD® AbSeq PCR1 primer	1
●	Library reverse primer 1–4	1 each
●	Library forward primer	1
●	Bead resuspension buffer	1
●	Elution buffer	1

BD® OMICS-One Dual Index Kit (-20 °C)		
Cap Color	Name	Quantity
●	Dual index forward primer 1–8	1 each
●	Dual index reverse primer 1–8	1 each

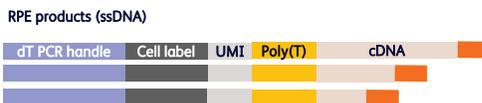
Workflows

WTA library amplification workflow

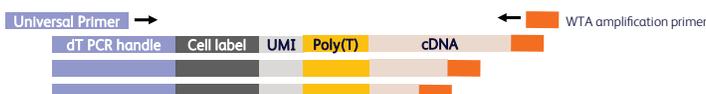
1.1 Random priming and extension (RPE) (page 13): Random priming on the bead.



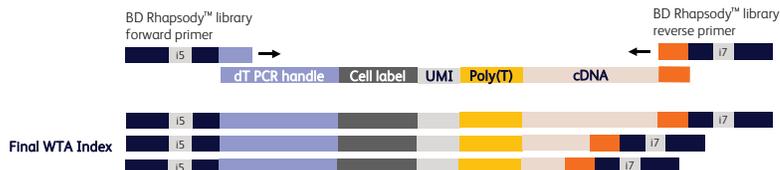
Denature off the RPE product.



1.4 RPE PCR (page 18): Amplify the RPE product.



1.6 WTA index PCR (page 22): Add adapters and indices.

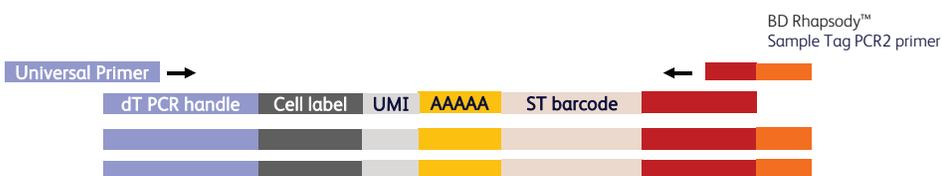


Sample Tag library amplification workflow

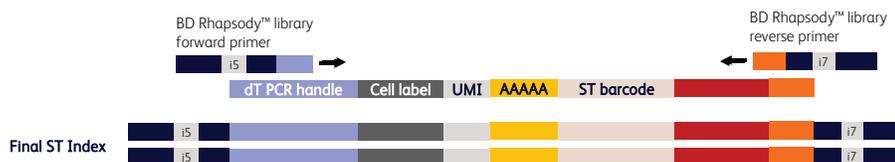
2.1 Sample Tag PCR1 (page 28): Sample Tag product is amplified in PCR1.



2.3 Sample Tag PCR2 (page 31): Sample Tag product is further amplified in PCR2.



2.5 Sample Tag index PCR (page 33): Add Sample Tag adapters and indices.



Required and recommended materials

Required reagents

Store the reagents at the storage temperature specified on the label.

Material	Supplier	Catalog no.
BD Rhapsody™ WTA Amplification Kit	BD Biosciences	633801
BD® OMICS-One Dual Index Kit [‡]	BD Biosciences	571899
Agencourt® AMPure® XP magnetic beads	Beckman Coulter	A63880
100% ethyl alcohol, molecular biology grade	Major supplier	–
Nuclease-free water	Major supplier	–

[‡] Recommended for unique dual indexing with high-throughput (more than eight) library preparation workflows.

Recommended consumables

Material	Supplier	Part number/Catalog no.
Pipettes (P10, P20, P200, P1000)	Major supplier	–
Low-retention, filtered pipette tips	Major supplier	–
0.2-mL PCR 8-strip tubes	Major supplier	–
Axygen® 96-Well PCR Microplates* Or, MicroAmp Optical 96-Reaction Plate*	Corning Thermo Fisher Scientific	PCR96HSC N8010560
MicroAmp Clear Adhesive Film*	Thermo Fisher Scientific	4306311
15-mL conical tube	Major supplier	–
DNA LoBind® tubes, 1.5 mL	Eppendorf	0030108051
Qubit™ Assay Tubes	Thermo Fisher Scientific	Q32856
Qubit™ dsDNA HS Assay Kit	Thermo Fisher Scientific	Q32851
Agilent High Sensitivity DNA Kit Or, Agilent High Sensitivity D1000 ScreenTape Agilent High Sensitivity D1000 Reagents Or, Agilent High Sensitivity D5000 ScreenTape Agilent High Sensitivity D5000 Reagents	Agilent Agilent Agilent Agilent Agilent	5067-4626 5067-5584 5067-5585 5067-5592 5067-5593

* Recommended for processing high-throughput (more than eight) library preparation workflows.

Equipment

Material	Supplier	Catalog no.
Microcentrifuge for 1.5–2.0-mL tubes	Major supplier	–
Microcentrifuge for 0.2-mL tubes	Major supplier	–
Vortexer	Major supplier	–
Digital timer	Major supplier	–
PCR thermal cycler	Major supplier	–
Eppendorf ThermoMixer® C*	Eppendorf	5382000023
6-tube magnetic separation rack for 1.5-mL tubes Or, 12-tube magnetic separation rack† Or, Invitrogen™ DynaMag™-2 magnet†	New England Biolabs New England Biolabs Thermo Fisher Scientific	S1506S S1509S 12321D
Low-profile magnetic separation stand for 0.2 mL, 8-strip tubes	V&P Scientific, Inc.	VP772F4-1
Magnetic Stand–96‡	Thermo Fisher Scientific	AM10027
Qubit™ 3.0 Fluorometer	Thermo Fisher Scientific	Q33216
Agilent® 2100 Bioanalyzer Or, Agilent® 4200 TapeStation System	Agilent Technologies Agilent Technologies	G2940CAG G2991AA
Heat block	Major supplier	–

* Two thermomixers are recommended. A heat block can be used for denaturation steps.

† Recommended for processing greater than six samples.

‡ Recommended for processing high-throughput (more than eight) library preparation workflows.

Best practices

Cell capture

- For best results, ensure that cells have high viability before proceeding with cell capture.

Bead handling

- When working with BD Rhapsody™ Enhanced Cell Capture Beads, use low-retention filtered tips and LoBind® tubes.



Never vortex the beads. Pipet-mix only.

- Store BD Rhapsody™ Enhanced Cell Capture Beads at 4 °C. Do not freeze.
- Bring Agencourt® AMPure® XP magnetic beads to room temperature (15–25 °C) before use. See the *AMPure® XP User's Guide* for information.

Libraries

- Sample Tag libraries can be sequenced separately or together with WTA mRNA libraries.

Master mix preparation

- Thaw reagents (except for enzymes) at room temperature.
- Keep enzymes at –25 °C to –15 °C until ready for use.
- Return reagents to correct storage temperature as soon as possible after preparing the master mix.

Supernatant handling

- Read this protocol carefully before beginning each section. Note which steps require you to keep supernatant to avoid accidentally discarding required products.
- Remove supernatants without disturbing AMPure® XP magnetic beads.
- Make and use fresh 80% ethyl alcohol within 24 hours. Adjust the volume of 80% ethyl alcohol depending on the number of libraries.

Additional documentation

- *BD Rhapsody™ HT Single-Cell Analysis System Single-Cell Capture and cDNA Synthesis Protocol* (doc ID 23-24252)
- *BD Rhapsody™ HT Xpress System Single-Cell Capture and cDNA Synthesis Protocol* (doc ID 23-24253)
- *BD Rhapsody™ System Single-Cell Labeling with BD® Single-Cell Multiplexing Kits Protocol* (doc ID 23-21340)
- *BD Rhapsody™ Sequence Analysis Pipeline User's Guide* (doc ID 23-24580)

Safety information

For safety information, refer to the *BD Rhapsody™ HT Single-Cell Analysis System Instrument User Guide* (doc ID 23-24607) or the *BD Rhapsody™ HT Xpress System Instrument User Guide for Scanner-Free Workflow* (doc ID 23-24256).

Time considerations

Station	Workflow	Timing	Stopping point and storage
WTA Library Amplification			
Pre-Amplification Workspace	Denature Sample Tags	10 minutes	4 °C up to 24 hours
Pre-Amplification Workspace	1.1 Random priming and extension (RPE) (page 13)	170 minutes	–
	1.2 RPE cleanup (page 16)		
Post-Amplification Workspace	1.4 RPE PCR (page 18)*	80 minutes	PCR overnight
	1.5 RPE PCR cleanup and quality check (page 19)		 <6 weeks at 4 °C or <6 months at –20 °C
	1.6 WTA index PCR (page 22)	65 minutes	PCR overnight
	1.7 WTA index PCR cleanup and quality check (page 23)		 <6 months at –20 °C
	(Optional) 1.8 Additional WTA index PCR cleanup (page 27)	25 minutes	 <6 months at –20 °C
* While the thermomixer program is running, proceed to the Sample Tag PCR1 task in the Sample Tag Library Amplification section.			
Sample Tag Library Amplification			
Post-Amplification Workspace	2.1 Sample Tag PCR1 (page 28)	65 minutes	PCR overnight
	2.2 Sample Tag PCR1 cleanup (page 30),		 <6 weeks at 4 °C or <6 months at –20 °C
	2.3 Sample Tag PCR2 (page 31)	60 minutes	PCR overnight
	2.4 Sample Tag PCR2 cleanup and quantify (page 31)		 <6 weeks at 4 °C or <6 months at –20 °C
	2.5 Sample Tag index PCR (page 33)	65 minutes	PCR overnight
	2.6 Sample Tag index PCR cleanup and quality check (page 35)		 <6 months at –20 °C
b. After PCR program has started, continue with RPE in the WTA Library Amplification section.			

Procedure

Continue this procedure after staining the antibodies as described in the *BD Rhapsody™ System Single-Cell Labeling with BD® Single-Cell Multiplexing Kits Protocol* or the *BD Rhapsody™ System Single-Cell Labeling with BD® Flex Single-Cell Multiplexing Kits Protocol*.

Perform the experiment on the BD Rhapsody™ Single-Cell Analysis system using either of the following guides for cell capture, reverse transcription, and Exonuclease treatment:

- *BD Rhapsody™ HT Single-Cell Analysis System Single-Cell Capture and cDNA Synthesis Protocol* (doc ID 23-24252)
- *BD Rhapsody™ HT Xpress System Single-Cell Capture and cDNA Synthesis Protocol* (doc ID 23-24253)

This protocol is intended for the whole transcriptome amplification library generation of cell inputs between 100 to 100,000 single cells, specifically resting PBMCs. For cell inputs between 100 and 5,000 single cells, follow the extra steps outlined in the additional cleanup sections.

Ensure that the intended total cell load is between 100 and 100,000 single cells for this protocol. Cell load below or above this recommendation might not be suitable for the current protocol configuration. Then proceed as described in the following procedure.

1. WTA library amplification

This procedure comprises the following tasks:

- [1.1 Random priming and extension \(RPE\) \(page 13\)](#)
- [1.2 RPE cleanup \(page 16\)](#)
- [1.3 Additional RPE cleanup for cell input <5,000 PBMC cells \(page 17\)](#)
- [1.4 RPE PCR \(page 18\)](#)
- [1.5 RPE PCR cleanup and quality check \(page 19\)](#)
- [1.6 WTA index PCR \(page 22\)](#)
- [1.7 WTA index PCR cleanup and quality check \(page 23\)](#)
- [1.8 Additional WTA index PCR cleanup \(page 27\)](#)

Before you begin

- Obtain Exonuclease I-treated and inactivated BD Rhapsody™ Enhanced Cell Capture Beads.
- Thaw reagents (except the enzymes) in the BD Rhapsody™ WTA Amplification Kit at room temperature (15–25 °C), then immediately place on ice.

1.1 Random priming and extension (RPE)

This section describes how to generate random priming products. First, random primers are hybridized to the cDNA on the BD Rhapsody™ Enhanced Cell Capture Beads, followed by extension with an enzyme.



Perform this procedure in the pre-amplification workspace. We recommend using a separate heat block for the 95 °C incubations.

1. Set a heat block to 95 °C and set two thermomixers to 37 °C and 25 °C, respectively.
2. In a new 1.5-mL LoBind® tube, pipet the following reagents.

Random primer mix

Cap	Component	For 1 library (µL)	For 1 library with 20% overage (µL)	For 4 libraries with 20% overage (µL)	For 8 libraries with 20% overage (µL)
●	WTA extension buffer	20.0	24.0	96.0	192.0
●	WTA extension primers	20.0	24.0	96.0	192.0
○	Nuclease-free water	134.0	160.8	643.2	1286.4
	Total	174.0	208.8	835.2	1670.4

3. Pipet-mix the random primer mix and keep at room temperature.
4. Briefly centrifuge the tube of Exonuclease I-treated BD Rhapsody™ Enhanced Cell Capture Beads, and then complete one of the following actions.
 - If you are using a subsample of the beads, proceed to the next step.
 - If you are using the entire sample of beads, skip to step 6.
5. (Optional) To subsample the Exonuclease I-treated BD Rhapsody™ Enhanced Cell Capture Beads:
 - a. Based on the expected number of viable cells captured on the beads in the final bead resuspension volume, determine the volume of beads to subsample for sequencing.
 - b. Completely resuspend the beads by pipet-mixing, then pipet the calculated volume of the bead suspension into a new 1.5-mL LoBind® tube.



The remaining Exonuclease I-treated beads can be stored in bead resuspension buffer at 4 °C for up to 1 year

6. Place the tube of Exonuclease I-treated beads in bead resuspension buffer on the 1.5-mL magnet for <2 minutes or until the supernatant is clear. Remove and discard the supernatant.
7. Remove the tube from the magnet and resuspend the beads in 75 µL of elution buffer. Pipet-mix 10 times to resuspend the beads. Label a new 1.5-mL tube as Sample Tag products.



If processing more than one library, we recommended performing the Sample Tag denaturation one library at a time.

8. Place the tube with beads in a 95 °C heat block for 5 minutes (no shaking).

9. Briefly centrifuge the tube, then place the tube on 1.5-mL magnet for <2 minutes or until the supernatant is clear.
 - a. Immediately remove the supernatant and transfer to the Sample Tag products tube from step 7.
 - b. To minimize Sample Tag contamination in the WTA library, ensure that all supernatant is removed from the tube.
 - c. Keep the supernatant tube on ice or at 4 °C for up to 24 hours until ready to proceed to [2.1 Sample Tag PCR1 \(page 28\)](#).
10. Remove the tube of BD Rhapsody™ Enhanced Cell Capture Beads from the magnet, and use a low-retention tip to pipet 200 µL of elution buffer into the tube. Pipet-mix 10 times to resuspend the beads.



If you are processing more than one library, beads in elution buffer can be stored on ice until all tubes had been denatured.

11. Briefly centrifuge the tube, then place the tube on a 1.5-mL magnet for <2 minutes or until the supernatant is clear. Remove and dispose of the supernatant.
12. Remove the tube with the BD Rhapsody™ Enhanced Cell Capture Beads from the magnet, and use a low-retention tip to pipet 87 µL of random primer mix into the tube. Pipet-mix 10 times to resuspend the beads.



Save the remaining volume of random primer mix for a second RPE. Keep random primer mix at room temperature.

13. Incubate the tube in the following order:
 - a. 95 °C in a heat block (no shaking) for 5 minutes.
 - b. Thermomixer at 1,200 rpm and at 37 °C for 5 minutes.
 - c. Thermomixer at 1,200 rpm and at 25 °C for 5 minutes.



Confirm “Time Mode” is set to “Time Control” before the program begins.

14. Briefly centrifuge the tube and keep it at room temperature.
15. Program the thermomixer.
 - a. 1,200 rpm and at 25 °C for 10 minutes.
 - b. 1,200 rpm and at 37 °C for 15 minutes.
 - c. 1,200 rpm and at 45 °C for 10 minutes.
 - d. 1,200 rpm and at 55 °C for 10 minutes.



Confirm “Time Mode” is set to “Time Control” before the program begins.

16. In a new 1.5-mL LoBind[®] tube, pipet the following reagents.

Extension enzyme mix

Cap	Component	For 1 library (μL)	For 1 library with 20% overage (μL)	For 4 libraries with 20% overage (μL)	For 8 libraries with 20% overage (μL)
	dNTP	8.0	9.6	38.4	76.8
	Bead RT/PCR enhancer	12.0	14.4	57.6	115.2
	WTA extension enzyme	6.0	7.2	28.8	57.6
	Total	26.0	31.2	124.8	249.6

17. Pipet 13 μL of the extension enzyme mix into the sample tube containing the beads (for a total volume of 100 μL) and keep at room temperature until ready.



Save the remaining volume of primer extension enzyme mix for a second RPE. Keep primer extension enzyme mix on ice.

18. Place the tube of extension enzyme mix with BD Rhapsody™ Enhanced Cell Capture Beads in the programmed thermomixer (see step 15).



While the thermomixer program is running, begin [2.1 Sample Tag PCR1 \(page 28\)](#).

19. Remove the tube after the program is finished. Place the tube in a 1.5-mL tube magnet for <2 minutes or until the supernatant is clear. Remove and discard the supernatant.

20. Remove the tube from the magnet and resuspend the beads in 205 μL of elution buffer using a P200 pipette.



If processing more than one library, we recommended performing the denaturation one library at a time.

21. To denature the random priming products off the beads, pipet to resuspend the beads.

- Incubate the sample at 95 °C in a heat block for 5 minutes (no shaking).
- Immediately after completing the 95 °C incubation, slightly open the lid of the tube to release air pressure within the tube.



Do not incubate for more than 5 minutes.

- Place the tube in a thermomixer at any temperature for 10 seconds at 1,200 rpm to resuspend the beads.

22. Place the tube in a 1.5-mL tube magnet for <2 minutes or until the supernatant is clear.
 - a. Immediately transfer 200 μ L of the supernatant containing the RPE product to a new 1.5-mL LoBind® tube.
 - b. Store supernatant containing the RPE product on ice.



If you are processing more than one library, supernatant containing RPE product can be stored on ice until all tubes had been denatured.

23. Repeat steps 12 through 21 to perform a second RPE.
24. Place the tube in a 1.5-mL tube magnet for <2 minutes or until the supernatant is clear.
 - a. Immediately transfer 200 μ L of the supernatant containing the RPE Product to the 1.5-mL LoBind® tube containing the supernatant from the first round of RPE (step 22) for a total of 400 μ L.
 - b. Store supernatant containing RPE product on ice. Discard the beads.
25. Immediately proceed to [1.2 RPE cleanup \(page 16\)](#).

1.2 RPE cleanup

This section describes how to perform a single-sided AMPure® cleanup, which removes primer dimers and other small molecular weight byproducts. The final product is purified single-stranded DNA.



Perform the purification in the pre-amplification workspace. Bring Agencourt® AMPure® XP magnetic beads to room temperature.

1. In a new 15-mL conical tube, prepare 10 mL of fresh 80% (v/v) ethyl alcohol by pipetting 8 mL of absolute ethyl alcohol to 2 mL of nuclease-free water. Vortex the tube for 10 seconds.



Make fresh 80% ethyl alcohol and use within 24 hours. The 80% ethyl alcohol volume should be adjusted depending on the number of libraries.

2. Vortex the AMPure® XP magnetic beads at high speed for 1 minute until the beads are fully resuspended.
3. If RPE product volume is <400 μ L, bring volume to 400 μ L with elution buffer. Pipet 720 μ L of AMPure® beads into the tube containing the 400 μ L of RPE product supernatant. Pipet-mix at least 10 times, then briefly centrifuge.
4. Incubate the suspension at room temperature for 10 minutes.
5. Place the tube on the 1.5-mL tube magnet for 5 minutes or until the supernatant is clear. Remove and discard the supernatant.
6. Keeping the tube on the magnet, gently pipet 1 mL of fresh 80% ethyl alcohol into the tube.
7. Incubate the sample on the magnet for 30 seconds. Remove and discard the supernatant.
8. Repeat steps 6 and 7 for a total of two ethyl alcohol washes.

9. Keeping the tube on the magnet, use a P20 pipette to remove and discard any residual supernatant from the tube.
 - a. For best results, briefly centrifuge the AMPure[®] beads while still wet and place the tube back on the magnet.
 - b. Remove and discard any excess ethanol that might collect at the bottom.
10. Air dry the beads at room temperature until no longer glossy (~15–20 minutes).

 Do not overdry the AMPure[®] beads after the ethanol washes. Overdried beads appear cracked.
11. Remove the tube from the magnet and resuspend the bead pellet in 40 µL of elution buffer. Pipet-mix the suspension at least 10 times until the beads are fully suspended.
12. Incubate the sample at room temperature for 2 minutes. Briefly centrifuge the tube to collect the contents at the bottom.
13. Place the tube on the magnet until the solution is clear, usually ~30 seconds.
14. Pipet the eluate (~40 µL) to a new PCR tube. This is the purified RPE product.
15. (Optional) For samples with low cell input (for example, starting with fewer than 5,000 PBMCs cell capture), proceed to [1.3 Additional RPE cleanup for cell input <5,000 PBMC cells \(page 17\)](#) for an additional round of AMPure[®] XP magnetic bead purification.
16. Keep on ice until ready to proceed with [1.4 RPE PCR \(page 18\)](#).



Additional RPE cleanup is only necessary when starting with <5,000 cells captured in the 8-lane cartridge. It is not necessary when processing <5,000 cells from subsampled beads.

1.3 Additional RPE cleanup for cell input <5,000 PBMC cells

1. To the tube from [1.2 RPE cleanup \(page 16\)](#), bring the purified RPE product volume up to 100 µL with nuclease-free water and transfer to a 1.5-mL LoBind[®] tube.

 The final volume must be exactly 100 µL to achieve the appropriate size selection of the purified RPE product.
2. Pipet-mix 10 times, then briefly centrifuge.
3. Pipet 180 µL of AMPure[®] beads into the tube containing 100 µL of eluted RPE product from the first round of purification.
4. Pipet-mix 10 times, then briefly centrifuge.
5. Incubate the suspension at room temperature for 5 minutes.
6. Place the suspension on the 1.5-mL tube magnet for 5 minutes or until the supernatant is clear. Remove and discard the supernatant.
7. Keeping the tube on the magnet, gently pipet 500 µL of fresh 80% ethyl alcohol into the tube.
8. Incubate the sample on the magnet for 30 seconds. Remove and discard the supernatant.
9. Repeat steps 7 and 8 for a total of two ethyl alcohol washes.

10. Keeping the tube on the magnet, use a small-volume pipette to remove and discard any residual supernatant from the tube.
11. Air dry the beads at room temperature until no longer glossy.
12. Remove the tube from the magnet and resuspend the bead pellet in 40 μL of elution buffer. Pipet-mix the suspension at least 10 times until the beads are fully suspended.
13. Incubate the sample at room temperature for 2 minutes. Briefly centrifuge the tube to collect the contents at the bottom.
14. Place the tube on the magnet until the solution is clear, usually ~30 seconds.
15. Pipet the eluate (~40 μL) to a new PCR tube. This is the purified RPE product after two rounds of purification.



Do not overdry the AMPure[®] beads after the ethanol washes. Overdried beads appear cracked.



STOPPING POINT: Store the RPE product in a LoBind[®] tube on ice or at 4 °C for up to 24 hours until [1.4 RPE PCR \(page 18\)](#).

1.4 RPE PCR

This section describes how to generate more RPE product through PCR amplification, resulting in multiple copies of each random-primed molecule.

1. In the pre-amplification workspace, in a new 1.5-mL LoBind[®] tube, pipet the following components.

RPE PCR mix

Cap	Component	For 1 library (μL)	For 1 library with 20% overage (μL)	For 4 libraries with 20% overage (μL)	For 8 libraries with 20% overage (μL)
○	PCR master mix	60	72	288	576
○	Universal oligo	10	12	48	96
○	WTA amplification primer	10	12	48	96
	Total	80	96	384	768

2. Pipet-mix the RPE PCR mix and keep on ice.
3. Add 80 μL of the RPE PCR mix to the tube with the 40 μL of purified RPE product. Pipet-mix 10 times to create the RPE PCR reaction mix.
4. Split the mix into two 0.2 mL PCR tubes with 60 μL of RPE PCR reaction mix per tube.

5. Bring the RPE PCR reaction mix to the post-amplification workspace and run the following PCR program.

RPE PCR program

Step	Cycles	Temperature	Time
Hot start	1	95 °C	3 min
Denaturation	Refer to the Recommended number of PCR cycles given in the following table*	95 °C	30 seconds
Annealing		60 °C	1 min
Extension		72 °C	1 min
Final extension	1	72 °C	2 min
Hold	1	4 °C	∞

* Recommended PCR cycles might require optimization for different cell types and cell number.

Recommended number of PCR cycles

Number of cells in RPE	Recommended PCR cycles for resting PBMCs
100	16
1,000–9,999	13
10,000	12
20,000	11
40,000	10
80,000–100,000	9

6. When the RPE PCR reaction is complete, briefly centrifuge to collect the contents at the bottom of the tubes.



STOPPING POINT: The PCR can run overnight.

1.5 RPE PCR cleanup and quality check

This section describes how to perform a single-sided AMPure® cleanup to remove unwanted small molecular weight products from the RPE products. The final product is purified double-stranded DNA (~200–2,000 bp).



Perform the purification in the post-amplification workspace.

1. Bring AMPure® XP magnetic beads to room temperature.
2. In a new 15-mL conical tube, prepare 5 mL of fresh 80% (v/v) ethyl alcohol by pipetting 4 mL of absolute ethyl alcohol to 1 mL of nuclease-free water. Vortex the tube for 10 seconds.



Make fresh 80% ethyl alcohol and use within 24 hours. The 80% ethyl alcohol volume should be adjusted depending on the number of libraries.

3. Vortex the AMPure® XP magnetic beads at high speed for 1 minute until the beads are fully resuspended.

4. Briefly centrifuge the tube with the RPE PCR product, then combine the two 60- μ L RPE PCR reactions into a new 1.5-mL tube.



The final volume must be exactly 120 μ L to achieve the appropriate size selection of the purified RPE PCR product. If the volume is less than 120 μ L, use elution buffer to achieve the final volume.

5. Pipet 120 μ L of AMPure[®] XP magnetic beads into the tube containing 120 μ L of RPE PCR product. Pipet-mix at least 10 times, then briefly centrifuge the samples.



Avoid getting AMPure[®] beads on the lid of the tube. Residual AMPure[®] beads and PCR mix buffer can negatively impact downstream results.

6. Incubate at room temperature for 5 minutes.
7. Place the 1.5-mL LoBind[®] tube on the magnet for 5 minutes or until the supernatant is clear. Remove and discard the supernatant.
8. Keeping the tube on the magnet, gently pipet 500 μ L of fresh 80% ethyl alcohol into the tube and incubate for 30 seconds. Remove and discard the supernatant without disturbing the beads.
9. Repeat step 8 once for a total of two ethyl alcohol washes.
10. Keeping the tube on the magnet, use a small-volume pipette to remove and discard any residual supernatant from the tube.
11. Air-dry the beads at room temperature for 5 minutes or until the beads no longer look glossy.



Do not overdry the AMPure[®] beads after the ethanol washes. Overdried beads appear cracked.

12. Remove the tube from the magnet and pipet 40 μ L of elution buffer into the tube to resuspend the bead pellet. Pipet-mix the suspension at least 10 times until the beads are fully suspended.
13. Incubate the sample at room temperature for 2 minutes. Briefly centrifuge the tube to collect the contents at the bottom.
14. Place the tube on the magnet until the solution is clear, usually ~30 seconds.
15. Pipet the eluate (~40 μ L) into a new 1.5-mL LoBind[®] tube. The RPE PCR product is ready for [1.6 WTA index PCR \(page 22\)](#).



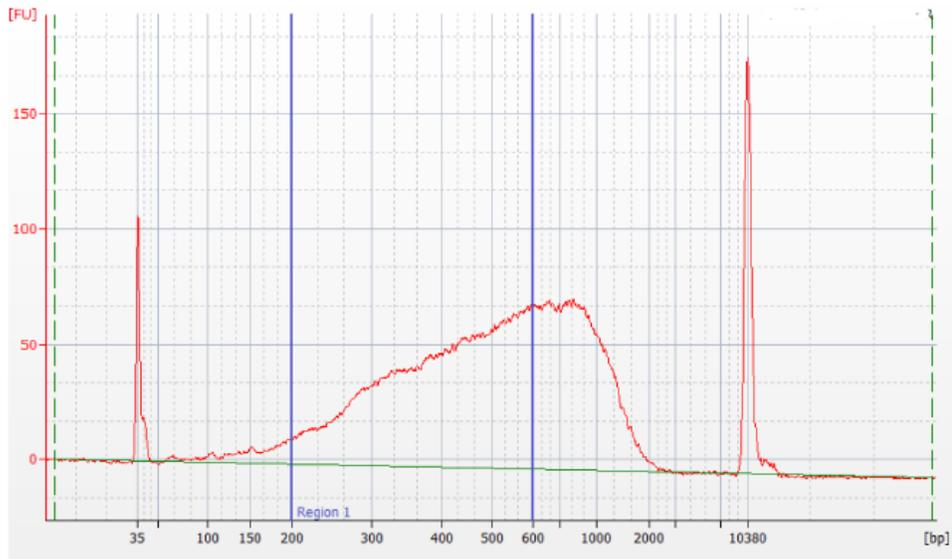
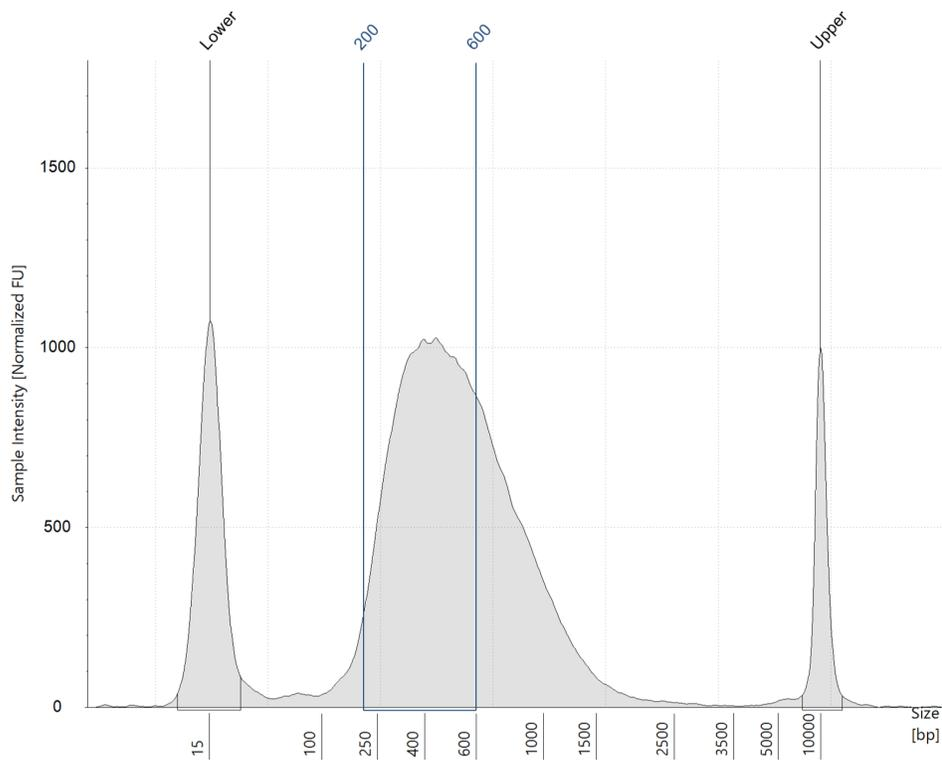
STOPPING POINT: The RPE PCR libraries can be stored at -20 °C for up to 6 months or 4 °C for up to 6 weeks.

16. Perform quality control of the RPE PCR products with the Agilent 2100 Bioanalyzer using the Agilent High Sensitivity DNA Kit or the Agilent 4200 TapeStation system using the Agilent High Sensitivity D1000 or D5000 ScreenTape assay.
 - a. The Bioanalyzer or TapeStation trace should show a broad peak from ~200 to 2,000 bp. Use the concentration from 200 to 600 bp to calculate how much template to add into Index PCR. Refer to the blue-boxed regions in the sample trace images in the following figures.

The Bioanalyzer or TapeStation is used to calculate molarity for the WTA library because of the distribution in fragment sizes for this library type.



Although there are products <200 bp (Sample Tag contamination product) and >600 bp, these products should be removed in the double-sided cleanup after the index PCR.

Figure 1 Sample bioanalyzer high-sensitivity DNA trace - RPE PCR product trace**Figure 2** Sample TapeStation high-sensitivity D5000 trace

1.6 WTA index PCR

This section describes how to generate mRNA libraries compatible with various sequencing platforms, by adding full-length sequencing adapters and indices through PCR.



Perform this procedure in the post-amplification workspace.



If additional unique or combinatorial indexing is needed, use the BD® OMICS-One Dual Index Kit primers.

1. Dilute the RPE PCR products from [1.5 RPE PCR cleanup and quality check \(page 19\)](#) with nuclease-free water until the concentration of the 200–600 bp peak is 2 nM. If the product concentration is <2 nM, do not dilute. Continue to the next step.

Example: If the Bioanalyzer measurement of the 200–600 bp peak is 6 nM, then dilute the sample threefold with nuclease-free water to 2 nM.

2. In a new 1.5-mL tube, pipet the following components to create the WTA index PCR mix.

WTA index PCR mix

Cap	Component	For 1 library (μL)	For 1 library with 20% overage (μL)	For 4 libraries with 20% overage (μL)	For 8 libraries with 20% overage (μL)
○	PCR master mix	25	30	120	240
●	Library forward primer	5	6	24	48
●	Library reverse primer 1–4*	5	6	–	–
○	Nuclease-free water	5	6	24	48
	Total	40	48	168	336

* For more than one library, use different library reverse primers for each library.

For sequencing on Illumina systems, refer to the Illumina guidelines for preparing libraries with balanced index combinations.

3. Pipet-mix the WTA index PCR mix and keep on ice.
4. In a new 0.2-mL PCR tube, combine WTA index PCR mix with diluted RPE PCR products as follows:
 - For one library: Combine 40 μL of WTA index PCR mix with 10 μL of 2 nM of RPE PCR product.
 - For multiple libraries: In separate tubes for each library, combine 35 μL of WTA Index PCR mix with 5 μL of the corresponding library reverse primer and 10 μL of 2 nM of RPE PCR products.
5. Pipet-mix 10 times.

6. Run the following PCR program.

Index PCR conditions for WTA

Step	Cycles	Temperature	Time
Hot start	1	95 °C	3 min
Denaturation	Refer to the Recommended number of PCR cycles shown in the following table*	95 °C	30 seconds
Annealing		60 °C	30 seconds
Extension		72 °C	30 seconds
Final extension	1	72 °C	1 min
Hold	1	4 °C	∞

* Cycle number varies based on the concentration of the RPE PCR product.

Recommended number of PCR cycles

Concentration of diluted RPE PCR products	Recommended number of PCR cycles
1 to <2 nM	9
2 nM	8

If the concentrations of diluted RPE PCR products are <1 nM, additional PCR cycles might be needed.



STOPPING POINT: The PCR can run overnight.

7. When the WTA index PCR is complete, briefly centrifuge to collect the contents at the bottom of the tubes.

1.7 WTA index PCR cleanup and quality check

This section describes how to perform a double-sided AMPure® cleanup for sequencing. The final product is purified double-stranded DNA with full-length adapter sequences.



Perform the purification in the post-amplification workspace.

1. Bring AMPure® XP magnetic beads to room temperature.
2. In a new 15-mL conical tube, prepare 5 mL of fresh 80% (v/v) ethyl alcohol by pipetting 4 mL of absolute ethyl alcohol to 1 mL of nuclease-free water. Vortex the tube for 10 seconds.
3. Vortex the AMPure® XP magnetic beads at high speed for 1 minute until the beads are fully resuspended.
4. Add 60 µL of nuclease-free water to the WTA Index PCR product for a final volume of 110 µL.
5. Transfer 100 µL of WTA Index PCR product into a new 0.2-mL PCR tube.



The volume must be exactly 100 µL.

6. Add 60 μL of AMPure[®] XP magnetic beads to the 0.2-mL PCR tube from the previous step. Pipet-mix at least 10 times, then briefly centrifuge.
7. Incubate the suspension at room temperature for 5 minutes.
8. Place the suspension on the 0.2-mL strip tube magnet for 3 minutes or until the supernatant is clear.
9. While the strip tube is still on the magnet, carefully remove and transfer the 160 μL of supernatant into a new 0.2 mL strip tube without disturbing the beads.
10. Pipet 15 μL of AMPure[®] beads into the 0.2 mL strip tube with supernatant (from step 9). Pipet-mix at least 10 times, then briefly centrifuge.



Discard the tubes with the pelleted AMPure[®] beads that contain long fragments.

11. Incubate the suspension at room temperature for 5 minutes.
12. Place the suspension on a 0.2-mL tube magnet for 3 minutes or until the supernatant is clear. Remove and discard the supernatant.
13. Keeping the tube on the magnet, gently pipet 200 μL of fresh 80% ethyl alcohol into the tube and incubate for 30 seconds. Discard the supernatant.
14. Repeat step 13 for a total of two ethyl alcohol washes.
15. Keeping the tube on the magnet, use a small-volume pipette to remove any residual supernatant from the tube.
16. Air-dry the beads at room temperature for 30 seconds.



Do not overdry the AMPure[®] beads after the ethanol washes. Overdried beads appear cracked.

17. Remove the tube from the magnet and pipet 30 μL of elution buffer into the tube. Pipet-mix at least 10 times to completely resuspend the AMPure[®] XP magnetic beads.
18. Incubate the sample at room temperature for 2 minutes.
19. Briefly centrifuge the tubes to collect the contents at the bottom.
20. Place the tube on the magnet until the solution is clear, usually ~30 seconds.
21. Pipet the eluate (~30 μL) into a new 1.5-mL LoBind[®] tube. The WTA Index PCR eluate is the final sequencing libraries.



STOPPING POINT: The Index PCR libraries can be stored at $-20\text{ }^{\circ}\text{C}$ for up to 6 months until sequencing.

22. Quantify and perform quality control of the Index PCR libraries with a Qubit™ Fluorometer using the Qubit™ dsDNA HS Assay and one of the following systems:
- The Agilent 2100 Bioanalyzer using the Agilent High Sensitivity DNA Kit
 - The Agilent 4200 TapeStation system using the Agilent High Sensitivity D1000 or D5000 ScreenTape assay



The expected concentration from the Qubit™ Fluorometer is >1 ng/ μ L. The Bioanalyzer trace should show a peak from ~ 250 to 1,000 bp. Refer to the sample trace images in the following figures.

Figure 3 Sample bioanalyzer high-sensitivity DNA trace–WTA index PCR product

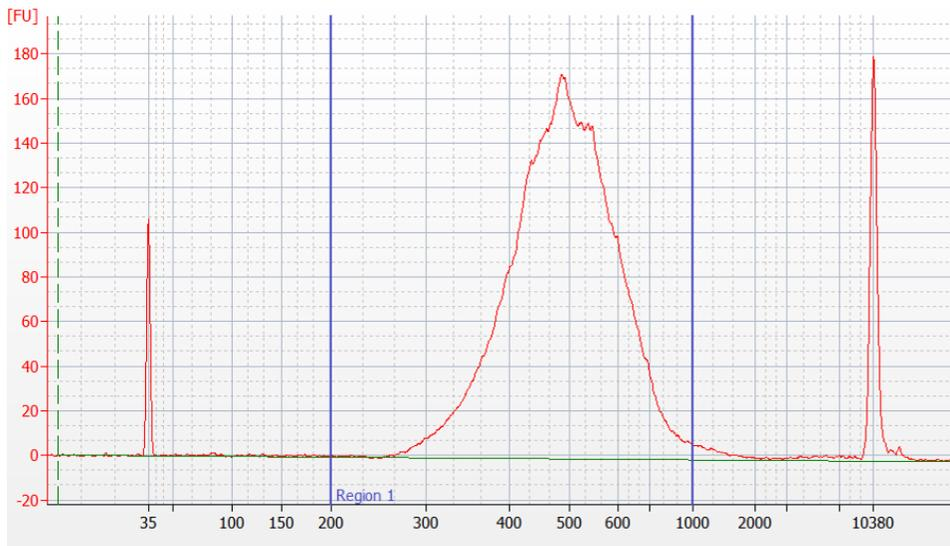
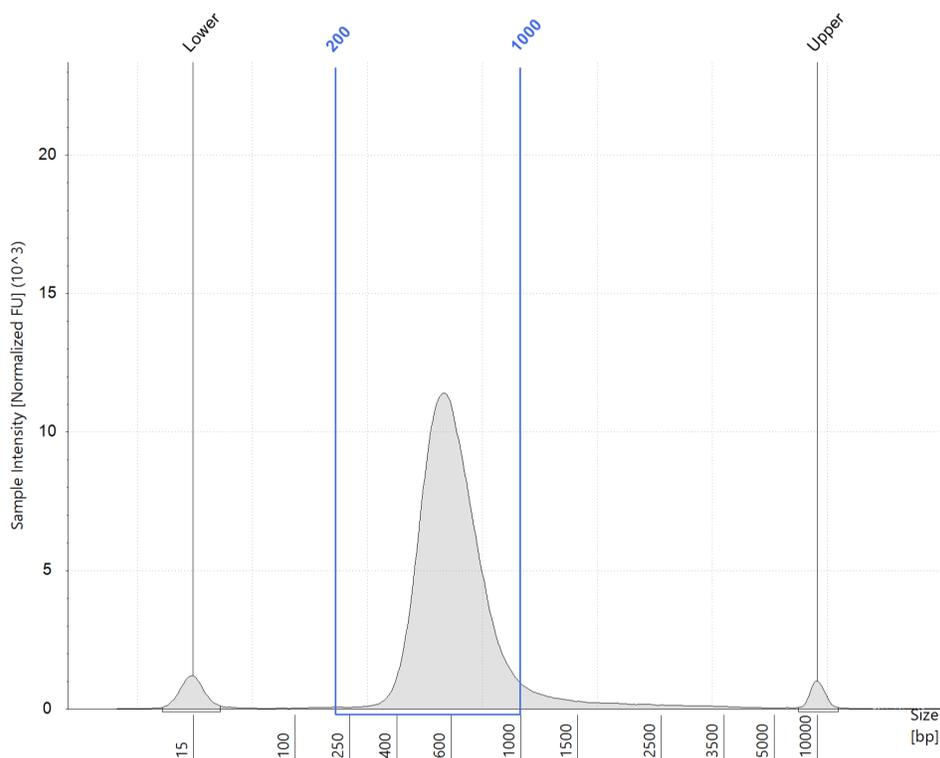
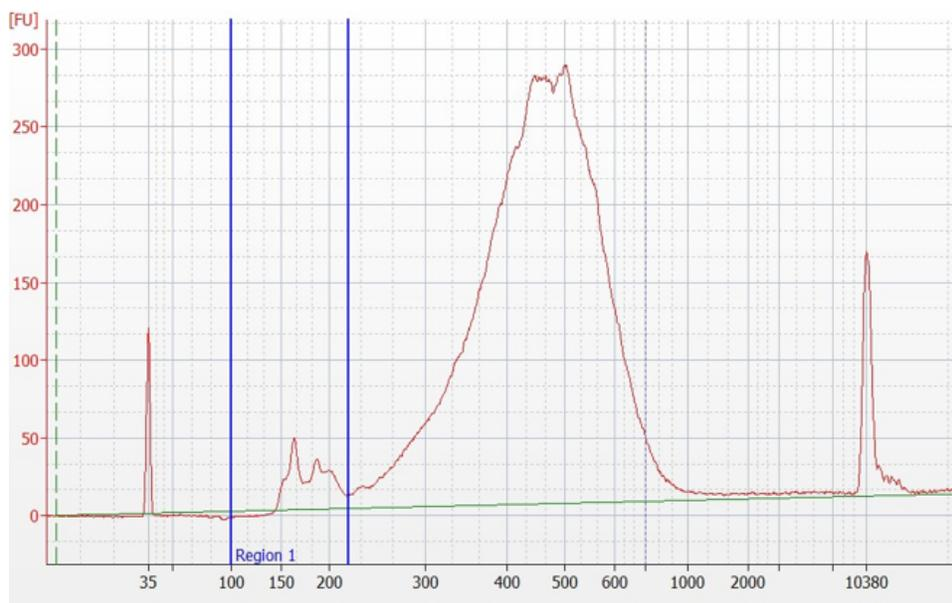


Figure 4 Sample TapeStation high-sensitivity D5000 trace–WTA index PCR product**Figure 5** Sample bioanalyzer high-sensitivity DNA trace for a WTA index PCR product with an observable <270 bp peak

If smaller products (~165 bp or ~270 bp) are observed (such as the peak shown in the preceding figure), we recommend a second round of AMPure® XP magnetic bead purification. See [1.8 Additional WTA index PCR cleanup \(page 27\)](#) for more information.

1.8 Additional WTA index PCR cleanup

If peaks <270 bp are observed, (as shown in the example [figure traces in 1.7 WTA index PCR cleanup and quality check \(page 25\)](#)), perform a second round of AMPure® XP magnetic bead purification.

1. To the tube from [step 21 in 1.7 WTA index PCR cleanup and quality check \(page 24\)](#), bring the total purified WTA index PCR eluate volume up to 100 µL with nuclease-free water. Pipet-mix 10 times, then briefly centrifuge.



The final volume must be exactly 100 µL to achieve the appropriate size selection of the WTA Index PCR product.

2. Pipet 75 µL of AMPure® beads into the tube containing 100 µL of eluted WTA index product from the first round of purification. Pipet-mix 10 times, then briefly centrifuge.
3. Incubate the suspension at room temperature for 5 minutes.
4. Place the suspension on the 1.5-mL tube magnet for 3 minutes or until the supernatant is clear.
5. Keeping the tube on the magnet, gently pipet 200 µL of fresh 80% ethyl alcohol into the tube and incubate the sample on the magnet for 30 seconds. Remove and discard the supernatant.
6. Repeat step 5 for a total of two ethyl alcohol washes.
7. Keeping the tubes on the magnet, use a small-volume pipette to remove and discard any residual supernatant from the tube.
8. Air-dry the beads at room temperature for 30 seconds.



Do not overdry the AMPure® beads after the ethanol washes. Overdried beads appear cracked.

9. Remove the tube from the magnet and resuspend the bead pellet in 30 µL of elution buffer. Pipet-mix the suspension at least 10 times until the beads are fully suspended.
10. Incubate the sample at room temperature for 2 minutes. Briefly centrifuge the tube to collect the contents at the bottom.
11. Place the tube on the magnet until the solution is clear, usually ~30 seconds.
12. Pipet the eluate (~30 µL) into a new 1.5-mL LoBind® tube. The WTA index PCR eluate is the final sequencing libraries.
13. Repeat the [last step in WTA index PCR cleanup and quality check \(page 25\)](#) to perform quality control.



STOPPING POINT: The Index PCR libraries can be stored at –20 °C for up to 6 months until sequencing.

2. Sample Tag library amplification

This section comprises the following tasks:

- [2.1 Sample Tag PCR1 \(page 28\)](#)
- [2.2 Sample Tag PCR1 cleanup \(page 30\)](#)
- [2.3 Sample Tag PCR2 \(page 31\)](#)
- [2.4 Sample Tag PCR2 cleanup and quantify \(page 31\)](#)
- [2.5 Sample Tag index PCR \(page 33\)](#)
- [2.6 Sample Tag index PCR cleanup and quality check \(page 35\)](#)

2.1 Sample Tag PCR1

This section describes how to amplify Sample Tag products through PCR.

1. In the pre-amplification workspace, pipet the following reagents into a new 1.5-mL LoBind® tube on ice.

Sample Tag PCR1 mix

Cap	Component	For 1 library (μL)	For 1 library with 20% overage (μL)	For 4 libraries with 20% overage (μL)	For 8 libraries with 20% overage (μL)
○	PCR master mix	100.0	120.0	480.0	960.0
○	Universal oligo	10.0	12.0	48.0	96.0
●	Sample Tag PCR1 primer	1.0	1.2	4.8	9.6
○	Nuclease-free water	21.0	25.2	100.8	201.6
	Total	132.0	158.4	633.6	1,267.2

2. Pipet-mix and keep on ice.
3. In a new 1.5-mL tube, pipet 132 μL of the Sample Tag PCR1 mix. Add 68 μL of the Sample Tag product from [step 9 in 1.1 Random priming and extension \(RPE\) \(page 14\)](#). Pipet-mix 10 times to create the Sample Tag PCR1 reaction mix. Do not vortex.
4. Pipet 50 μL Sample Tag PCR1 reaction mix into each of four 0.2-mL PCR tubes. Transfer any residual mix to one of the tubes.
5. Bring the Sample Tag PCR1 reaction mix to the post-amplification workspace.
6. Run the following PCR program on the thermal cycler.

PCR1 conditions for Sample Tag panel

Step	Cycles	Temperature	Time
Hot start	1	95 °C	3 min
Denaturation	8–16*	95 °C	30 seconds
Annealing		60 °C	30 seconds
Extension		72 °C	1 min
Final extension	1	72 °C	5 min
Hold	1	4 °C	∞

* Recommended PCR cycles might need to be optimized for different cell types and cell number.

Recommended number of PCR cycles

Number of cells in PCR1	Recommended PCR cycles for resting PBMCs
100	16
1,000	14
2,500	13
5,000	12
10,000	11
20,000	10
40,000	9
80,000–100,000	8



STOPPING POINT: The PCR can be run overnight.

- After PCR has started, continue with [1.1 Random priming and extension \(RPE\) \(page 13\)](#).
- After PCR, briefly centrifuge the tubes.
- Pipet-mix and combine the four reactions into a new 1.5-mL LoBind[®] tube labeled Sample Tag PCR1. Keep the tube on ice and proceed to [2.2 Sample Tag PCR1 cleanup \(page 30\)](#).

2.2 Sample Tag PCR1 cleanup

This section describes how to perform a single-sided AMPure® cleanup to remove primer dimers from the Sample Tag PCR1 products. The final product is purified double-stranded DNA.



Perform the purification in the post-amplification workspace.

1. Bring AMPure® XP magnetic beads to room temperature.
2. In a new 15-mL conical tube, prepare 5 mL of fresh 80% (v/v) ethyl alcohol by combining 4 mL absolute ethyl alcohol, molecular biology grade, with 1 mL nuclease-free water. Vortex the tube for 10 seconds to mix.



Make fresh 80% ethyl alcohol and use it within 24 hours. The 80% ethyl alcohol volume should be adjusted depending on the number of libraries.

3. Vortex on high speed for 1 minute until the AMPure® beads are fully resuspended.
4. Briefly centrifuge the Sample Tag PCR1 products.



The final volume must be exactly 200 μ L to achieve the appropriate size selection of the purified Sample Tag PCR1 product. If the volume is less than 200 μ L, use elution buffer to achieve the final volume.

5. To 200 μ L of Sample Tag PCR1 products, pipet 360 μ L AMPure® beads (from [2.1 Sample Tag PCR1 \(page 28\)](#)).
6. Pipet-mix 10 times. Incubate at room temperature for 5 minutes.
7. Place the suspension on the 1.5-mL tube magnet for 5 minutes or until the supernatant is clear. Remove and discard the supernatant.
8. Keeping the tube on the magnet, gently pipet 500 μ L of fresh 80% ethyl alcohol into the tube and incubate for 30 seconds. Remove and discard the supernatant.
9. Repeat step 8 once for a total of two ethyl alcohol washes.
10. Keeping the tube on the magnet, use a small-volume pipette to remove and discard any residual supernatant from the tube.
11. Air-dry the beads at room temperature for 5 minutes.



Do not overdry the AMPure® beads after the ethanol washes. Overdried beads appear cracked.

12. Remove the tube from the magnet and resuspend the bead pellet in 30 μ L of elution buffer. Pipet-mix the suspension at least 10 times until the beads are fully suspended. Small clumps do not affect performance.
13. Incubate at room temperature for 2 minutes and briefly centrifuge.
14. Place the tube on the magnet until the solution is clear, usually ~30 seconds.
15. Pipet the eluate (~30 μ L) into a new 1.5-mL LoBind® tube (purified Sample Tag PCR1 products).



STOPPING POINT: Store at 2–8 °C before proceeding within 24 hours or at –25 °C to –15 °C for up to 6 months

2.3 Sample Tag PCR2

This section describes how to amplify Sample Tag products through PCR. The PCR primers include partial sequencing adapters that enable the additions of full-length sequencing indices during index PCR.

1. In the pre-amplification workspace, pipet reagents into a new 1.5-mL LoBind[®] tube on ice.

Sample Tag PCR2 mix

Cap	Component	For 1 library (μL)	For 1 library with 20% overage (μL)	For 4 libraries with 20% overage (μL)	For 8 libraries with 20% overage (μL)
○	PCR master mix	25.0	30.0	120.0	240.0
○	Universal oligo	2.0	2.4	9.6	19.2
●	Sample Tag PCR2 primer	3.0	3.6	14.4	28.8
○	Nuclease-free water	15.0	18.0	72.0	144.0
	Total	45.0	54.0	216.0	432.0

2. Pipet-mix the Sample Tag PCR2 mix and keep on ice.
3. Pipet 45 μL of Sample Tag PCR2 mix into one 0.2-mL PCR tube for each library.
4. Bring the Sample Tag PCR2 mix to the post-amplification workspace.
5. Pipet 5 μL of PCR1 products into 45 μL Sample Tag PCR2 mix to create the Sample Tag PCR2 reaction mix. Pipet-mix 10 times and briefly centrifuge.
6. Run the following PCR program on the thermal cycler.

PCR2 conditions for Sample Tag

Step	Cycles	Temperature	Time
Hot start	1	95 °C	3 min
Denaturation	10*	95 °C	30 seconds
Annealing		66 °C	30 seconds
Extension		72 °C	1 min
Final extension	1	72 °C	5 min
Hold	1	4 °C	∞

* Recommended PCR cycles might need to be optimized for different cell types and cell number.



STOPPING POINT: The PCR can run overnight.

2.4 Sample Tag PCR2 cleanup and quantify

This section describes how to perform a single-sided AMPure[®] cleanup to remove primer dimers from the Sample Tag PCR2 products. The final product is purified double-stranded DNA.



Perform PCR2 purification in the post-amplification workspace.

1. Bring AMPure® XP magnetic beads to room temperature.
2. In a new 5.0-mL LoBind® tube, prepare 5 mL fresh 80% (v/v) ethyl alcohol by combining 4 mL absolute ethyl alcohol, molecular biology grade, with 1 mL of nuclease-free water. Vortex the tube for 10 seconds to mix.



Make fresh 80% ethyl alcohol and use it within 24 hours. The 80% ethyl alcohol volume should be adjusted depending on the number of libraries.

3. Vortex AMPure® XP beads at high speed for 1 minute until beads are fully resuspended.
4. Briefly centrifuge the Sample Tag PCR2 products.



The final volume must be exactly 50 µL to achieve the appropriate size selection of the purified Sample Tag PCR2 product. If the volume is less than 50 µL, use elution buffer to achieve the final volume.

5. Pipet 60 µL of AMPure® XP beads into 50.0 µL of Sample Tag PCR2 product.
6. Pipet-mix 10 times and incubate at room temperature for 5 minutes.
7. Place the tube on the strip tube magnet for 3 minutes or until the supernatant is clear. Remove and discard the supernatant.
8. Keeping the tube on the magnet, gently pipet 200 µL of fresh 80% ethyl alcohol into the tube and incubate for 30 seconds. Remove and discard the supernatant.
9. Repeat step 8 for a total of two ethyl alcohol washes.
10. Keeping the tube on the magnet, use a small-volume pipette to remove and discard any residual supernatant from the tube.
11. Air-dry the beads at room temperature for 2 minutes.



Do not overdry the AMPure® beads after the ethanol washes. Overdried beads appear cracked.

12. Remove the tube from the magnet and resuspend the bead pellet in 30 µL of elution buffer. Pipet-mix the suspension at least 10 times until the beads are fully resuspended.
13. Incubate at room temperature for 2 minutes and briefly centrifuge to collect the contents at the bottom.
14. Place the tube on the magnet until the solution is clear, usually ~30 seconds.
15. Pipet the eluate (~30 µL) into a new 1.5-mL LoBind® tube. The Sample Tag PCR2 product is ready for Sample Tag index PCR.



STOPPING POINT: Store at 2–8 °C before proceeding on the same day, or at –25 °C to –15 °C for up to 6 months.

16. Estimate the concentration by quantifying 2 µL of the final sequencing library with a Qubit™ Fluorometer using the Qubit™ dsDNA HS Assay Kit. Follow the manufacturer's instructions.
17. Dilute an aliquot of the products with nuclease-free water to 0.1–1.1 ng/µL.

2.5 Sample Tag index PCR

This section describes how to generate Sample Tag libraries compatible with various sequencing platforms, by adding full-length sequencing adapters and indices through PCR.



For cell capture samples from multiple lanes, the same reverse primer can be used to label all the library types from one lane (for example, WTA and Sample Tag from Lane 1 can both be labeled with reverse primer 1, while WTA and Sample Tag from Lane 2 can be labeled with reverse primer 2, and so on). The kit provides 4 indexing primers, and can label all sample combinations from up to 4 lanes for the same sequencing run.



If additional unique or combinatorial indexing is needed, use the BD[®] OMICS-One Dual Index Kit primers.

1. In the pre-amplification workspace, pipet reagents into a new 1.5-mL LoBind[®] tube on ice.

Sample Tag index PCR mix

Cap	Component	For 1 library (μL)	For 1 library with 20% overage (μL)	For 4 libraries with 20% overage (μL)	For 8 libraries with 20% overage (μL)
○	PCR master mix	25.0	30.0	120.0	240.0
●	Library forward primer	2.0	2.4	9.6	19.2
●	Library reverse primer 1–4*	2.0	2.4	–	–
○	Nuclease-free water	18.0	21.6	86.4	172.8
	Total	47.0	56.4	216.0	432.0

* For more than one library, use different library reverse primers for each library.

For sequencing on Illumina systems, refer to the Illumina guidelines for preparing libraries with balanced index combinations.

2. Pipet-mix the Sample Tag index PCR mix, and keep on ice.
3. Bring the Sample Tag index PCR mix to post-amplification workspace.
4. In a new 0.2 mL PCR tube, combine Sample Tag index PCR mix with diluted Sample Tag PCR2 products as follows based on the number of libraries:
 - For one Sample Tag library, pipet 3.0 μL of 0.1–1.1 ng/μL Sample Tag PCR2 product into 47.0 μL Sample Tag index PCR mix (from [2.4 Sample Tag PCR2 cleanup and quantify \(page 31\)](#)). See the Concentration Index PCR input for Sample Tag libraries (ng/μL) in [Recommended number of PCR cycles \(page 34\)](#).
 - For multiple libraries, combine 45.0 μL of Sample Tag index PCR mix with 2 μL of corresponding library reverse primer, and then pipet 3.0 μL of 0.1–1.1 ng/μL Sample Tag PCR2 product into the PCR tube.
5. Gently vortex, and briefly centrifuge.

6. Run the following PCR program on the thermal cycler.

Index PCR conditions for Sample Tag

Step	Cycles	Temperature	Time
Hot start	1	95 °C	3 min
Denaturation	Refer to the Recommended number of PCR cycles in the following table*	95 °C	30 seconds
Annealing		60 °C	30 seconds
Extension		72 °C	30 seconds
Final extension	1	72 °C	1 min
Hold	1	4 °C	∞

* Cycle number varies based on the concentration of the Sample Tag PCR2 product.

Recommended number of PCR cycles

Concentration Index PCR input for Sample Tag libraries (ng/μL)	Recommended number of PCR cycles
0.5–1.1	6
0.25–0.5	7
0.1–0.25	8



STOPPING POINT: The PCR can run overnight.

2.6 Sample Tag index PCR cleanup and quality check

This section describes how to perform a single-sided AMPure[®] cleanup to remove primer dimers from the Sample Tag Index PCR products. The final product is purified double-stranded DNA with full-length adapter sequences.



Perform Index PCR purification in the post-amplification workspace.

1. Bring AMPure[®] XP magnetic beads to room temperature.
2. In a new 15-mL LoBind[®] tube, prepare 5 mL fresh 80% (v/v) ethyl alcohol by combining 4 mL absolute ethyl alcohol, molecular biology grade, with 1 mL of nuclease-free water. Vortex the tube for 10 seconds to mix.



Make fresh 80% ethyl alcohol, and use it within 24 hours. The 80% ethyl alcohol volume should be adjusted depending on the number of libraries.

3. Vortex the AMPure[®] XP magnetic beads at high speed for 1 minute until the beads are fully resuspended.
4. Briefly centrifuge all the index PCR products.



The final volume must be exactly 50 μ L to achieve the appropriate size selection of the purified Sample Tag index PCR products. If the volume is less than 50 μ L, use elution buffer to achieve the final volume.

5. Pipet 40 μ L of AMPure[®] beads into the 0.2 mL PCR tube containing 50.0 μ L of the Sample Tag index PCR product.
6. Pipet-mix 10 times and incubate at room temperature for 5 minutes.
7. Place the tube on the strip tube magnet for 3 minutes or until the supernatant is clear. Remove and discard the supernatant.
8. Keeping the tube on the magnet, gently add 200 μ L of fresh 80% ethyl alcohol into the tube, and incubate the sample on the magnet for 30 seconds. Discard the supernatant.
9. Repeat step 8 for a total of two ethyl alcohol washes.
10. Keeping the tube on the magnet, use a small-volume pipette to remove and discard the residual supernatant from the tube.
11. Air-dry the beads at room temperature for 2 minutes.



Do not overdry the AMPure[®] beads after the ethanol washes. Overdried beads appear cracked.

12. Remove the tube from the magnet and resuspend the bead pellet in 30 μ L of elution buffer. Pipet-mix the suspension at least 10 times until the beads are fully suspended.
13. Incubate at room temperature for 2 minutes. Briefly centrifuge the tube to collect the contents at the bottom.
14. Place the tube on the magnet until the solution is clear, usually ~30 seconds.

15. Pipet the entire eluate (~30 μ L) into a new 1.5-mL LoBind® tube. The index PCR product is ready for sequencing.



STOPPING POINT: Store at $-25\text{ }^{\circ}\text{C}$ to $-15\text{ }^{\circ}\text{C}$ for up to 6 months until sequencing.

16. Perform quality control before freezing samples.

- a. Estimate the concentration by quantifying with a Qubit™ Fluorometer using the Qubit™ dsDNA HS Kit to obtain an approximate concentration of PCR products to dilute for quantification on an Agilent 2100 Bioanalyzer or an Agilent 4200 TapeStation system using the Agilent High Sensitivity D1000 or D5000 ScreenTape Assay. Follow the manufacturer's instructions. The expected concentration of the libraries is $>1.5\text{ ng}/\mu\text{L}$.



The Sample Tag library should show a peak of ~276 bp. Exact size may vary due to instrument or sample purification efficiency. Select the region size between 200 and 500 bp.

Figure 6 Sample bioanalyzer high-sensitivity DNA trace—Sample Tag index PCR product

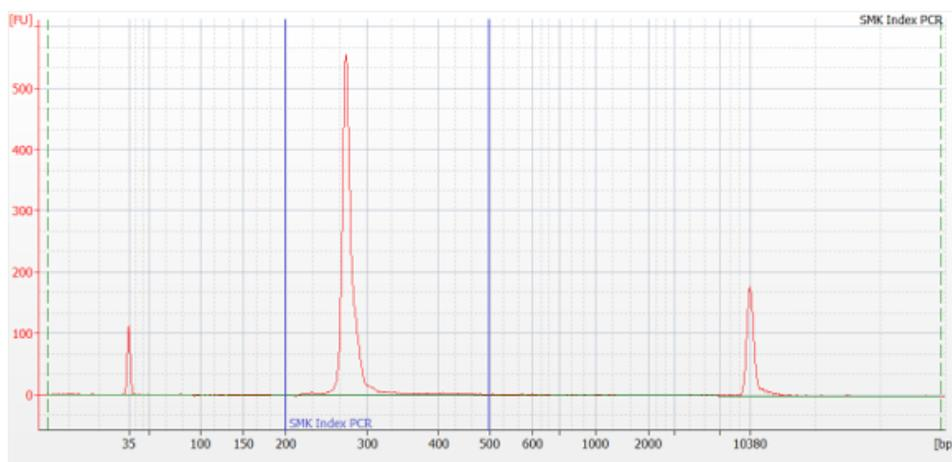
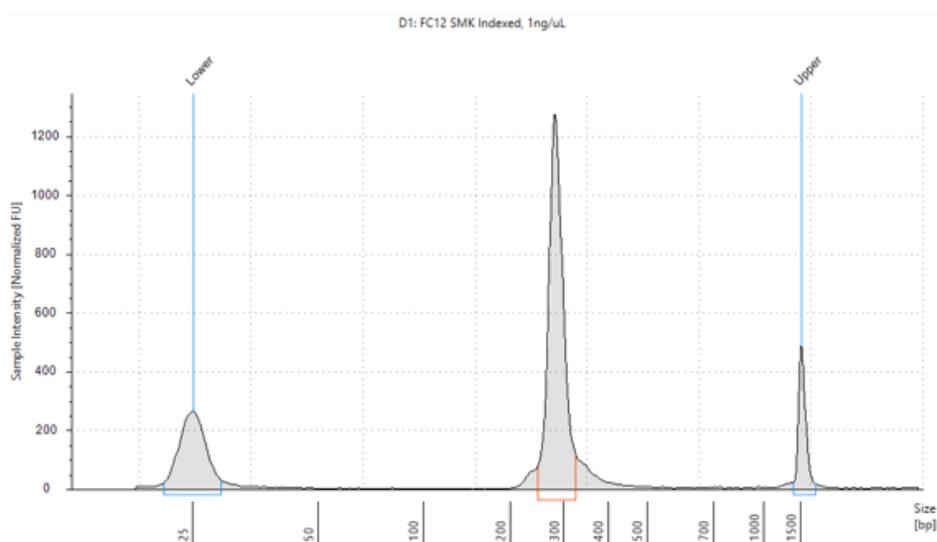


Figure 7 Sample TapeStation high-sensitivity D1000 trace—Sample Tag index PCR product



Sequencing

The sequencing depth for each library is dependent on application. For cell-type clustering, shallow sequencing is sufficient. However, for in-depth analysis such as comparison across multiple samples, deep sequencing is advised. We recommend meeting the requirement for recursive substitution error correction (RSEC) sequencing depth of ≥ 6 to reach the threshold of sequencing saturation where most molecules of the library have been recovered, approximately 80%. The RSEC sequencing depth and sequencing saturation are both reported by the analysis pipeline. The actual sequencing reads/cell required to achieve this depth can vary as it is dependent on the gene expression levels, number of cells, and sequencing run quality. The following reads/cell are recommended for WTA mRNA and Sample Tag libraries.

Read requirements for libraries

Gene panel	Read requirement for data analysis
BD Rhapsody™ WTA	20,000–100,000 reads/cell
Sample Tag	600 reads/cell

Required parameters

Parameter	Requirement
Platform	Illumina*
Paired-end reads	Recommend Read 1: 51 cycles; Read 2: 71 cycles; Index 1 (i5): 8 cycles; Index 2 (i7): 8 cycles
PhiX	1% recommended
Analysis	See the <i>BD® Single-Cell Multiomics Bioinformatics Handbook</i>

* To review Illumina Index 1 (i7) sequences, see the [Appendix \(page 42\)](#).

Sequencing recommendations

- For a NextSeq High or Mid Output run and MiniSeq High or Mid Output run, load the flow cell at a concentration between 1.5 and 1.8 pM with 1% PhiX for a sequencing run.
- For Novaseq:

Sequencing platform	Cycles	Recommended loading concentration
NovaSeq 6000 S Prime (Single Lane)	2×50, 2×100, 2×150, 2×250*	180–250 pM (XP workflow)
NovaSeq 6000 S Prime (Single Flow Cell)	2×50, 2×100, 2×150, 2×250*	350–650 pM (standard workflow)
NovaSeq 6000 S1 (Single Lane)	2×50, 2×100, 2×150*	180–250 pM (XP workflow)
NovaSeq 6000 S1 (Single Flow Cell)	2×50, 2×100, 2×150*	350–650 pM (standard workflow)
NovaSeq 6000 S2 (Single Flow Cell)	2×50, 2×100, 2×150*	350–650 pM (standard workflow)
NovaSeq 6000 S4 (Single Lane)	2×100, 2×150	180–250 pM (XP workflow)
NovaSeq 6000 S4 (Single Flow Cell)	2×100, 2×150	350–650 pM (standard workflow)
NovaSeq X 10B	2×100, 2×150	Contact local Field Application Specialist (FAS)

* NovaSeq 100 cycle kit (v1.0 or v1.5) can be used. The 100-cycle kit contains enough reagents for up to 130 cycles.

- For other sequencing platforms (e.g. Element AVITI System), follow the manufacturer's sequencing recommendations.

Sequencing depth can vary depending on whether the sample contains high- or low-content RNA cells. For resting PBMCs, we recommend:

- 10,000 reads per cell for shallow sequencing. Genes per cell and UMI per cell detected is generally lower, but this can be useful for cell type identification.
- 20,000–50,000 reads per cell for moderate sequencing.
- 100,000 reads per cell for highly saturated deep sequencing to identify the majority of UMIs in the library.



To determine the ratio of BD Rhapsody™ WTA mRNA library to Sample Tag library to pool for sequencing, use the sequencing calculator available by contacting your local Field Application Specialist (FAS) or scomix@bdscomix.bd.com.

Sequencing analysis pipeline

Contact customer support at scomix@bdscomix.bd.com for access to the latest whole transcriptome sequencing analysis pipeline.

Troubleshooting

Library preparation

Observation	Possible causes	Recommended solutions
PCR2 product yield too low.	PCR1 and PCR2 primers might have been swapped by mistake.	Ensure the correct primers are used for each step.
	cDNA synthesis might have failed due to incomplete washing of lysis buffer.	Avoid leaving behind lysis buffer or bubbles after removing lysis buffer from the tube during bead wash after retrieval from the cartridge. Use new tubes for each wash step, as described in the protocol.
	cDNA synthesis might have failed due to thermomixer not shaking during reverse transcription.	Samples need to be on the thermomixer in shake mode. Where applicable, ensure that a SmartBlock™ Thermoblock is installed on the thermomixer for 1.5-mL tubes so that the reaction can proceed at the designated temperature.
	Thermal cycler mis-programming.	Ensure that the correct thermal cycling program is used.
	Too few PCR1 cycles.	Optimize the number of PCR cycles for the specific sample type.
	Incorrect volume of AMPure® XP magnetic beads used during PCR2 cleanup.	Use the specified volume of AMPure® XP beads.
	Incorrect solution or incorrect concentration of 80% ethyl alcohol used for washing AMPure® XP magnetic beads, resulting in premature elution of PCR products from beads.	Use 80% ethyl alcohol for washing AMPure® XP beads.
Low yield of RPE-PCR.	Cell number lower than expected.	Repeat PCR using the RPE PCR product for additional cycles. Alternatively, increase index PCR cycles.
Index PCR BioAnalyzer trace of WTA library has 264 bp peak.	Sample Tag library contamination in mRNA library.	If peak takes up high percentage of sequencing reads (manifests as lower reads/cell than expected for WTA library, alongside higher reads/cell than expected for Sample Tag), perform a second round of AMPure® purification according to 1.8 Additional WTA index PCR cleanup (page 27) .
Low yield of indexing PCR.	Input DNA not high enough or cycle number too low.	Repeat indexing PCR with higher cycle number. Alternatively, if RPE-PCR product was diluted before adding to indexing PCR, repeat indexing PCR with less or no dilution.

Observation	Possible causes	Recommended solutions
Index PCR Bioanalyzer trace of WTA library shows large amount of product larger than 600 bp.	Over-amplification during indexing PCR.	Repeat indexing PCR with lower cycle number. Alternatively, repeat indexing with diluted RPE-PCR product.
Final sequencing product size too large	Over-amplification during index PCR or input amount of PCR2 products too high.	Repeat the index PCR with a lower input of PCR2 products.
	Upper and lower markers on the Agilent Bioanalyzer or TapeStation are incorrectly called.	Ensure that markers are correct. Follow manufacturer's instructions.
	Incorrect volume of Agencourt® AMPure® XP magnetic beads used.	Use volume specified in protocol.
Yield of Sample Tag library too low after index PCR (<1 ng/μL).	Sample Tag labeling incubation time too short.	Ensure that the cells were labeled with Sample Tags correctly and that the correct incubation time was used.
	PCR1 and PCR2 primers swapped.	Ensure that correct primer is used for each step.
	Only one primer (library forward or library reverse primer) added to index PCR mix.	Ensure that both the library forward primer and library reverse primer are added to the index PCR mix, and repeat index PCR.
	Too few index PCR cycles.	Increase the number of index PCR cycles.
Lower number of reads/cell than expected from mRNA.	264 bp or ~160 bp products taking reads from mRNA library.	If noise peak is seen in the 264 bp or ~160 bp range, perform a second round of AMPure® purification according to 1.8 Additional WTA index PCR cleanup (page 27) .
Expected size of Sample Tag products is too short (<276 bp).	Upper and lower markers on the Agilent Bioanalyzer or Agilent TapeStation are incorrectly called.	Ensure that the markers are correct. Follow the manufacturer's instructions.
	Inefficient Sample Tag labeling.	Ensure that the cells were labeled with Sample Tags correctly and that the correct incubation time was used.
	Sample Tags were not amplified in PCR steps due to incorrect primers used.	Perform PCR2 again. See 2.3 Sample Tag PCR2 (page 31) . Analyze products using the Agilent Bioanalyzer or the Agilent TapeStation and look for a ~182 bp peak that corresponds to Sample Tag PCR2 products.

Sequencing

Observation	Possible causes	Recommended solutions
Over-clustering on the Illumina flow cell due to under-estimation of the library.	Inaccurate measurement of the library concentration.	Quantitate library according to instructions in protocol.
Low sequencing quality.	Insufficient PhiX.	Use the recommended concentration of PhiX with the library to be sequenced.
	Suboptimal cluster density, or library denaturation, or both.	See troubleshooting in Illumina documentation.
High proportion of undetermined Sample Tag calls in sequencing results.	Insufficient sequencing of the Sample Tag Library.	<ol style="list-style-type: none"> 1. Set pooled samples of the same cell type to 120 reads/cell. 2. Set pooled samples of different cell types to 600 reads/cell. 3. Repeat sequencing of Sample Tag library. If issue persists, contact your local Field Application Specialist (FAS) or scomix@bdscomix.bd.com.
	Insufficient washes after labeling cells with Sample Tags.	Follow the washing steps in this protocol.
	BD Rhapsody™ Cartridge overloaded with cells.	Follow the cell loading steps in the <i>BD Rhapsody™ Single-Cell Analysis System Instrument User Guide</i> .

Appendix

Oligonucleotides in BD Rhapsody™ Whole Transcriptome Analysis Amplification Kit

The following table lists the sequences of all oligonucleotides included in the BD Rhapsody™ Whole Transcriptome Analysis Amplification Kit (Catalog No. 633801). Note that the BD® AbSeq primer and BD Rhapsody™ WTA amplification primer have the same sequence.

Oligonucleotide	Use	Part/ catalog no.	Sequence (5'–3')
BD Rhapsody™ Universal Oligo	Forward primer for WTA RPE PCR, Sample Tag PCR1 and PCR2, and BD® AbSeq PCR1	650000074	ACACGACGCTCTCCGATCT
BD Rhapsody™ Sample Tag PCR1 Primer	Reverse primer for Sample Tag PCR1	91-1088	GTTGTCAAGATGCTACCGTT
BD Rhapsody™ Sample Tag PCR2 Primer	Reverse primer for Sample Tag PCR2	91-1089	CAGACGTGTGCTCTCCGATCTGTTGTCAAGATGCTACCGTT
BD Rhapsody™ WTA Extension Primers	Random primers for WTA RPE	91-1115	TCAGACGTGTGCTCTCCGATCTNNNNNNNNN
BD Rhapsody™ WTA Amplification Primer	Reverse primer for WTA RPE PCR	91-1116	CAGACGTGTGCTCTCCGATCT
BD Rhapsody™ Library Forward Primer	Forward primer for WTA, Sample Tag, and BD® AbSeq Index PCR	91-1085	AATGATACGGCGACCACCGAGATCTACACTATAGCCT ACACTCTTCCCTACACGACGCTCTTCCGATC*T
BD Rhapsody™ Library Reverse Primer 1	WTA, Sample Tag, and BD® AbSeq Index PCR	650000080	CAAGCAGAAGACGGCATAACGAGATAGCGTAGCGTGA CTGGAGTTCAGACGTGTGCTCTTCCGATC*T
BD Rhapsody™ Library Reverse Primer 2		650000091	CAAGCAGAAGACGGCATAACGAGATCAGCCTCGGTGA CTGGAGTTCAGACGTGTGCTCTTCCGATC*T
BD Rhapsody™ Library Reverse Primer 3		650000092	CAAGCAGAAGACGGCATAACGAGATTGCCTCTTGTGA CTGGAGTTCAGACGTGTGCTCTTCCGATC*T
BD Rhapsody™ Library Reverse Primer 4		650000093	CAAGCAGAAGACGGCATAACGAGATTCCTCTACGTGA CTGGAGTTCAGACGTGTGCTCTTCCGATC*T

Forward index name	i5 bases for sample sheet NovaSeq, MiSeq, HiSeq 2000/2500	i5 bases for sample sheet iSeq, MiniSeq, NexSeq, HiSeq 3000/4000
BD Rhapsody™ Library Forward Primer	TATAGCCT	AGGCTATA

Reverse index name	i7 bases for sample sheet
BD Rhapsody™ Library Reverse Primer 1 (N709)	GCTACGCT
BD Rhapsody™ Library Reverse Primer 2 (N710)	CGAGGCTG
BD Rhapsody™ Library Reverse Primer 3 (N711)	AAGAGGCA
BD Rhapsody™ Library Reverse Primer 4 (N712)	GTAGAGGA

BD® OMICS-One Dual Index Kit sequences

Forward index name	i5 bases for sample sheet NovaSeq, MiSeq, HiSeq 2000/2500	i5 bases for sample sheet iSeq, MiniSeq, NexSeq, HiSeq 3000/4000
Dual index forward primer 1	TATAGCCT	AGGCTATA
Dual index forward primer 2	ATAGAGGC	GCCTCTAT
Dual index forward primer 3	CCTATCCT	AGGATAGG
Dual index forward primer 4	GGCTCTGA	TCAGAGCC
Dual index forward primer 5	AGGCGAAG	CTTCGCCT
Dual index forward primer 6	TAATCTTA	TAAGATTA
Dual Index Forward Primer 7	CAGGACGT	ACGTCCTG
Dual Index Forward Primer 8	GTA CTGAC	GTCAGTAC

Reverse index name	i7 bases for sample sheet
Dual index reverse primer 1	ATTACTCG
Dual index reverse primer 2	TCCGGAGA
Dual index reverse primer 3	CGCTCATT
Dual index reverse primer 4	GAGATTCC
Dual index reverse primer 5	ATTCAGAA
Dual index reverse primer 6	GAATTCGT
Dual index reverse primer 7	CTGAAGCT
Dual index reverse primer 8	TAATGCGC

Sample Tag sequences

Human Sample Tag sequences

Each Human Sample Tag is a human universal antibody conjugated with a unique oligonucleotide sequence to allow for sample identification. Each Sample Tag has common 5' and 3' ends, with the Sample Tag sequence between them:

- GTTGTCAAGATGCTACCGTTCAGAG(**Sample Tag sequence**)AAAAAAAAAAAAAAAAAAAAAAAAAAAA

Sample Tag	Sample Tag sequence
Sample Tag 1–Human	ATTCAAGGGCAGCCGCGTCACGATTGGATACGACTGTTGGACCGG
Sample Tag 2–Human	TGGATGGGATAAGTGCCTGATGGACCGAAGGGACCTCGTGGCCGG
Sample Tag 3–Human	CGGCTCGTGCTGCGTCTCAAGTCCAGAACTCCGTGTATCCT
Sample Tag 4–Human	ATTGGGAGGCTTTCGTACCGCTGCCGCCACCAGGTGATACCCGCT
Sample Tag 5–Human	CTCCCTGGTGTCAATACCCGATGTGGTGGGCAGAATGTGGCTGG
Sample Tag 6–Human	TTACCCGCAGGAAGACGTATAACCCCTCGTGCCAGGCGACCAATGC
Sample Tag 7–Human	TGTCTACGTCGGACCGCAAGAAGTGAGTCAGAGGCTGCACGCTGT
Sample Tag 8–Human	CCCCACCAGTTGCTTTGTCGGACGAGCCCGCACAGCGCTAGGAT
Sample Tag 9–Human	GTGATCCGCGCAGGCACACATACCGACTCAGATGGGTTGTCCAGG
Sample Tag 10–Human	GCAGCCGGCGTCGTACGAGGCACAGCGGAGACTAGATGAGGCCCC
Sample Tag 11–Human	CGCGTCCAATTTCCGAAGCCCCGCCCTAGGAGTCCCTGCGTGC
Sample Tag 12–Human	GCCCATTCAATGCACCCGCCAGTGATCGACCCTAGTGGAGCTAAG

Mouse Immune Sample Tag sequences

Each Mouse Immune Sample Tag is an Anti-Mouse CD45, Clone 30-F11 antibody conjugated with a unique oligonucleotide sequence to allow for sample identification. Each Sample Tag has common 5' and 3' ends, with the Sample Tag sequence between them:

- GTTGTCAAGATGCTACCGTTCAGAG(**Sample Tag sequence**)AAAAAAAAAAAAAAAAAAAAAAAAAAAA

Sample Tag	Sample Tag sequence
Sample Tag 1–Mouse Immune	AAGAGTCGACTGCCATGTCCCCTCCGCGGGTCCGTGCCCCCAAG
Sample Tag 2–Mouse Immune	ACCGATTAGGTGCGAGGCGCTATAGTCGTACGTCGTTGCCGTGCC
Sample Tag 3–Mouse Immune	AGGAGGCCCGCGTGAGAGTGATCAATCCAGGATACATTCCCGTC
Sample Tag 4–Mouse Immune	TTAACCGAGGCGTGAGTTTGGAGCGTACCGGCTTTGCGCAGGGCT
Sample Tag 5–Mouse Immune	GGCAAGGTGTCACATTGGGCTACCGCGGGAGGTCGACCAGATCCT
Sample Tag 6–Mouse Immune	GCGGGCACAGCGGCTAGGGTGTTCCGGGTGGACCATGGTTCAGGC
Sample Tag 7–Mouse Immune	ACCGGAGGCGTGTGTACGTGCGTTTCGAATCCTGTAAGCCCACC
Sample Tag 8–Mouse Immune	TCGCTGCCGTGCTTCATTGTCGCCGTTCTAACCTCCGATGTCTCG
Sample Tag 9–Mouse Immune	GCCTACCCGCTATGCTCGTCGGCTGGTTAGAGTTTACTGCACGCC
Sample Tag 10–Mouse Immune	TCCCATTGGAATCACGAGGCCGGGTGCGTTCTCCTATGCAATCCC
Sample Tag 11–Mouse Immune	GGTTGGCTCAGAGGCCCGAGGCTGCGGACGTCGTCGGACTCGCGT
Sample Tag 12–Mouse Immune	CTGGGTGCCTGGTCGGGTTACGTCGGCCCTCGGGTCGCGAAGGTC

BD® Flex SMK sequences

Each Flex Sample Tag is an anti-PE antibody conjugated with a unique oligonucleotide sequence to allow for sample identification. Each Sample Tag has common 5' and 3' ends, with the Sample Tag sequence between them:

- GTTGTCAAGATGCTACCGTTCAGAG(**Sample Tag sequence**)AAAAAAAAAAAAAAAAAAAAAAAAAAAA

BD® Flex Single-Cell Multiplexing Kit A (Catalog No. 633849)



Not for use together with Hu SMK Tags 1–6 (Catalog No. 633781) in the same experiment.

Sample Tag	Sample Tag sequence	Notes
Sample Tag 1–Flex	ATTCAAGGGCAGCCGCGTCACGATTGGATACGACTGTTGGACCGG	Barcode sequence is the same as human SMK Sample Tag 1
Sample Tag 2–Flex	TGGATGGGATAAGTGC GTGATGGACCGAAGGGACCTCGTGGCCGG	Barcode sequence is the same as human SMK Sample Tag 2
Sample Tag 3–Flex	CGGCTCGTGCTGCGTCGTCTCAAGTCCAGAACTCCGTGTATCCT	Barcode sequence is the same as human SMK Sample Tag 3
Sample Tag 4–Flex	ATTGGGAGGCTTTCGTACCGCTGCCGCCACCAGGTGATACCCGCT	Barcode sequence is the same as human SMK Sample Tag 4
Sample Tag 5–Flex	CTCCCTGGTGTTCAATACCCGATGTGGTGGGCAGAATGTGGCTGG	Barcode sequence is the same as human SMK Sample Tag 5
Sample Tag 6–Flex	TTACCCGCAGGAAGACGTATACCCCTCGTGCCAGGCGACCAATGC	Barcode sequence is the same as human SMK Sample Tag 6

BD[®] Flex Single-Cell Multiplexing Kit B (Catalog No. 633850)

Not for use together with Hu SMK Tags 7–12 (Catalog No. 633781) in the same experiment.

Sample Tag	Sample Tag sequence	Notes
Sample Tag 7–Flex	TGTCTACGTCCGACCGCAAGAAGTGAGTCAGAGGCTGCACGCTGT	Barcode sequence is the same as human SMK Sample Tag 7
Sample Tag 8–Flex	CCCCACCAGGTTGCTTTGTCGGACGAGCCCGCACAGCGCTAGGAT	Barcode sequence is the same as human SMK Sample Tag 8
Sample Tag 9–Flex	GTGATCCGCGCAGGCACACATACCGACTCAGATGGGTTGTCCAGG	Barcode sequence is the same as human SMK Sample Tag 9
Sample Tag 10–Flex	GCAGCCGGCGTCGTACGAGGCACAGCGGAGACTAGATGAGGCCCC	Barcode sequence is the same as human SMK Sample Tag 10
Sample Tag 11–Flex	CGCGTCCAATTTCCGAAGCCCCGCCCTAGGAGTCCCCTGCGTGC	Barcode sequence is the same as human SMK Sample Tag 11
Sample Tag 12–Flex	GCCCATTCATTGCACCCGCCAGTGATCGACCCTAGTGGAGCTAAG	Barcode sequence is the same as human SMK Sample Tag 12

BD® Flex Single-Cell Multiplexing Kit C (Catalog No. 633851)



Not for use together with Ms SMK Tags 1–6 (Catalog No. 633793) in the same experiment.

Sample Tag	Sample Tag sequence	Notes
Sample Tag 13–Flex	AAGAGTCGACTGCCATGTCCCCTCCGCGGGTCCGTGCCCCCAAG	Barcode sequence is the same as mouse SMK Sample Tag 1
Sample Tag 14–Flex	ACCGATTAGGTGCGAGGCGCTATAGTCGTACGTCGTTGCCGTGCC	Barcode sequence is the same as mouse SMK Sample Tag 2
Sample Tag 15–Flex	AGGAGGCCCGCGTGAGAGTGATCAATCCAGGATACATTCCCGTC	Barcode sequence is the same as mouse SMK Sample Tag 3
Sample Tag 16–Flex	TTAACCGAGGCGTGAGTTTGGAGCGTACCGGCTTTGCGCAGGGCT	Barcode sequence is the same as mouse SMK Sample Tag 4
Sample Tag 17–Flex	GGCAAGGTGTCACATTGGGCTACCGCGGGAGGTGCACCAGATCCT	Barcode sequence is the same as mouse SMK Sample Tag 5
Sample Tag 18–Flex	GCGGGCACAGCGGCTAGGGTGTCCGGGTGGACCATGGTTCAGGC	Barcode sequence is the same as mouse SMK Sample Tag 6

BD[®] Flex Single-Cell Multiplexing Kit D (Catalog No. 633852)

Not for use together with Ms SMK Tags 7–12 (Catalog No. 633793) in the same experiment.

Sample Tag	Sample Tag sequence	Notes
Sample Tag 19–Flex	ACCGGAGGCGTGTGTACGTGCGTTTCGAATCCTGTAAGCCCACC	Barcode sequence is the same as mouse SMK Sample Tag 7
Sample Tag 20–Flex	TCGCTGCCGTGCTTCATTGTGCGCCGTCTAACCTCCGATGTCTCG	Barcode sequence is the same as mouse SMK Sample Tag 8
Sample Tag 21–Flex	GCCTACCCGCTATGCTCGTCGGCTGGTTAGAGTTTACTGCACGCC	Barcode sequence is the same as mouse SMK Sample Tag 9
Sample Tag 22–Flex	TCCCATTCGAATCACGAGGCCGGGTGCGTTCTCCTATGCAATCCC	Barcode sequence is the same as mouse SMK Sample Tag 10
Sample Tag 23–Flex	GGTTGGCTCAGAGGCCCCAGGCTGCGGACGTCGCGGACTCGCGT	Barcode sequence is the same as mouse SMK Sample Tag 11
Sample Tag 24–Flex	CTGGGTGCCTGGTCGGGTTACGTGCGCCCTCGGGTCGCGAAGGTC	Barcode sequence is the same as mouse SMK Sample Tag 12

Contact Information

Becton, Dickinson and Company

BD Biosciences

155 North McCarthy Boulevard

Milpitas, California 95035 USA

bdbiosciences.com

scomix@bdscomix.bd.com