

 BD Rhapsody™ System
Extended-Lysis Single-Cell Capture and
cDNA Synthesis with
BD Rhapsody™ Express Single-Cell Analysis
System
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](https://www.bd.com/patents).

BD, the BD Logo, BD Rhapsody and Pharmingen are trademarks of Becton, Dickinson and Company or its affiliates. All other trademarks are the property of their respective owners. © 2025 BD. All rights reserved.

Regulatory information

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

History

Revision	Date	Change made
23-24981(01)	2025-09	Initial release.

Contents

Introduction	4
Symbols	4
Required and recommended materials	4
Required reagents	4
Recommended consumables	5
Equipment	6
Preparation list	7
Best practices	9
Before you begin	9
Priming and treating the BD Rhapsody™ Cartridge	10
Counting and preparing a single-cell suspension for cartridge loading	11
Loading cells in the cartridge	13
Preparing BD Rhapsody™ Enhanced Cell Capture Beads	14
Loading and washing BD Rhapsody™ Enhanced Cell Capture Beads	15
Lysing cells	16
Retrieving BD Rhapsody™ Enhanced Cell Capture Beads	17
Washing BD Rhapsody™ Enhanced Cell Capture Beads	18
Performing reverse transcription	19
Treating the sample with Exonuclease I	20
Troubleshooting	21

Introduction

This protocol describes cell loading in the BD Rhapsody™ Cartridge and single-cell capture with the BD Rhapsody™ Express Single-Cell Analysis System.

For complete instrument procedures, safety information, and workflow, see the *BD Rhapsody™ Express Single-Cell Analysis System Instrument User Guide*.

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.
	Biological hazard.

Required and recommended materials

For a complete list of materials, see the instrument user guide.

Required reagents

Material	Supplier	Catalog no.
BD Rhapsody™ Enhanced Cartridge Reagent Kit V3	BD Biosciences	667052
BD Rhapsody™ Cartridge Kit	BD Biosciences	633733
BD Rhapsody™ cDNA Kit	BD Biosciences	633773
Absolute ethyl alcohol, molecule biology grade	Major supplier	–
Nuclease-free water	Major supplier	–

Recommended consumables

Material	Supplier	Catalog no.
Gilson™ PIPETMAN™ Tipack™ Filter Tips, 100-1200 µL for BD Rhapsody™ P1200M pipette	Thermo Fisher Scientific	F171803G
Gilson™ PIPETMAN™ Tipack™ Filter Tips, 500-5000 µL for BD Rhapsody™ P5000M pipette	Thermo Fisher Scientific	F161370G
Falcon® Tube with Cell Strainer Cap	Corning	352235
DNA LoBind® Tubes, 1.5-mL	Eppendorf	30108051
DNA LoBind® Tubes, 5.0-mL ^a	Eppendorf	30108310
Low-retention, filtered pipette tips (10 µL, 200 µL, 1000 µL)	Major supplier	–
Pre-moistened cleaning wipes with 70% ethyl alcohol or 70% isopropyl alcohol.	Major supplier	–
Lint-free wipes	Major supplier	–
a. These are the Bead Retrieval Tubes to be used with the BD Rhapsody™ Express instrument.		

Equipment

Material	Supplier	Catalog no.
BD Rhapsody™ Express Instrument ^a	BD Biosciences	633702
BD Rhapsody™ P1200M pipette ^a	BD Biosciences	633704
BD Rhapsody™ P5000M pipette ^a	BD Biosciences	633705
Cell Counter	Major supplier	–
Large magnetic separation stand	V&P Scientific, Inc.	VP 772FB-1
Clear acrylic cylinder adapter for 15-mL tube adapter ^b	V&P Scientific, Inc.	VP 772FB-1A
Microcentrifuge for 1.5–2.0-mL tubes	Major supplier	–
Centrifuge and rotor with adapters for 5mL Falcon® tubes and 15-mL tubes	Major supplier	–
Eppendorf ThermoMixer® C	Eppendorf	5382000023
SmartBlock™ Thermoblock 1.5-mL	Eppendorf	5360000038
Plate Shaker for cartridge workflow: SmartBlock™ plates (for ThermoMixer C) Or, Eppendorf MixMate® Or, MicroPlate Genie™	Eppendorf Eppendorf Scientific Industries, Inc.	5363000039 22674200 SI-0400
Water bath OR incubator at 37 °C	Major supplier	–
Pipettes (P10, P20, P200, P1000)	Major supplier	–
Vortexer	Major supplier	–
Digital timer	Major supplier	–
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
a. Part of the BD Rhapsody™ Single-Cell Analysis system. Items can be ordered separately. b. Holds 5-mL LoBind tube in magnet.		

Preparation list

Item		BD Part Number	Preparation and Handling	Storage
Equilibrate to room temperature				
Cartridge Wash Buffer 1		650000060	Equilibrate to room temperature 30 minutes before setting up reverse transcription (RT). Centrifuge briefly. Keep on ice until ready.	4 °C
Cartridge Wash Buffer 2		650000061		-20 °C
	RT Buffer	650000067		
	dNTP	650000077		
	RT 0.1 M DTT	650000068		
	Nuclease-Free Water	650000076		
	10X Exonuclease I Buffer	650000071		
	Bead Resuspension Buffer	650000066		
Place on ice				
BD Rhapsody™ Enhanced Cell Capture Beads			Centrifuge briefly. Keep on ice until ready.	4 °C
Sample Buffer		650000062		
Lysis Buffer		650000064		
Bead Wash Buffer		650000065		
1.0 M DTT		650000063		
Leave at -25 °C to -15 °C				
	Bead RT/PCR Enhancer	91-1082	Centrifuge briefly before adding to mix.	-20 °C
	RNase Inhibitor	650000078		
	Reverse Transcriptase	700026321		
	Exonuclease I	650000072		
Obtain				
BD Rhapsody™ Cartridge Kit		633733	—	Ambient
2 mM Calcein AM			Thaw and protect from light.	-20 °C

Item	BD Part Number	Preparation and Handling	Storage
0.3 mM DRAQ7™		Protect from light.	4 °C
100% ethyl alcohol			
Ice bucket			
1.5 mL DNA LoBind® tubes			
INCYTO™ disposable hemocytometer			
Cell suspension			
Set up			
Thermomixer at 80 °C			
Thermomixer at 37 °C			

Best practices

- Always use low-retention filtered pipette tips and LoBind tubes.
- Perform single-cell capture and cDNA synthesis in a pre-amplification workspace.
- Prepare cells as close to cell loading as possible.
- Change pipette tips before every pipetting step.
- Keep reagents on ice unless instructed otherwise.
- To ensure an airtight seal with the BD Rhapsody™ P1200M and P5000M pipettes, hold the pipette with one hand and slightly twist the pipette to firmly seat a pipette tip on the pipette shaft.

Before you begin

- Visually inspect the Lysis Buffer for any precipitation. If precipitation is present, incubate the Lysis Buffer at room temperature (15–25 °C) for 1 hour. Invert to mix but do not vortex. Once the solution is clear, place the Lysis Buffer on ice.
- Open the DTT tube while holding the tube vertically. The solution is overlain with an inert/non-oxygen-containing gas, and a non-vertical tube will allow the inert gas to pour off. If not loading 4 or 8 lanes at the same time, after opening the DTT tube once, seal and store the tube at –20 °C or aliquot for single use in 0.5-mL microcentrifuge tubes and store at –20 °C.
- If cell preparation takes 4 hours or longer, begin preparing cells before cartridge preparation.
- For guidance on preparing a single-cell suspension, see *Preparing Single-Cell Suspensions Protocol*.
- If your biological sample contains red blood cell contamination, red blood cell lysis is required. See *Preparing Single-Cell Suspensions Protocol*.

Priming and treating the BD Rhapsody™ Cartridge

Prime and treat the BD Rhapsody™ Cartridge. For detailed instructions, see the instrument user guide.

Express instrument slider	Position
Front	Waste
Side	0

Step no.	Material to load	Volume (μL)	P1200M pipette mode	Incubation at room temperature
1	100% ethyl alcohol	700	Prime/Treat	—
2	Air	700	Prime/Treat	—
3	Room temp. Cartridge Wash Buffer 1	700	Prime/Treat	1 min
4	Air	700	Prime/Treat	—
5	Room temp. Cartridge Wash Buffer 1	700	Prime/Treat	10 min
6	Air	700	Prime/Treat	—
7	Room temp. Cartridge Wash Buffer 2	700	Prime/Treat	≤4 hr

Counting and preparing a single-cell suspension for cartridge loading

For detailed instructions, see the instrument user guide.

- 1 Treat the cells with viability stain(s), and count. Order of accurate counting:
 - Manual counting with fluorescence
 - Automated counting with fluorescence
 - Automated counting with Trypan Blue Stain and brightfield
 - Manual counting with Trypan Blue Stain and brightfield
- 2 Determine the desired number of cells to capture in the BD Rhapsody™ Cartridge . See the instrument user guide for a table containing estimated multiplet rates based on the number of captured cells on retrieved BD Rhapsody™ Enhanced Cell Capture Beads.
- 3 Determine the pooling ratio of samples to load onto the cartridge. For example, if two samples were labeled using the BD Rhapsody™ Single-Cell Multiplexing Kit, and the samples will be pooled in equal proportion, the pooling ratio for each sample is 0.5. If only one sample is used, the pooling ratio is 1.
- 4 Calculate the volume, V , for each sample needed to prepare the pooled single-cell suspension:

$$V = N \times P \times 1.36 / C$$

where:

V = volume of sample needed (μL)

N = desired number of captured cells in cartridge

P = pooling ratio

C = total cell concentration (cells/ μL)

Example

On a BD Rhapsody™ Cartridge, you want to capture 10,000 cells that are pooled equally of Sample A and Sample B.

N = desired number of captured cells in cartridge = 10,000

P_A = sample A pooling ratio = 0.5

P_B = sample B pooling ratio = 0.5

C_A = sample A total cell concentration = 200 cells/ μL

C_B = sample B total cell concentration = 400 cells/ μL

Volume of sample A needed = 10,000 cells $\times$ 0.5 $\times$ 1.36 / 200cells/ μL = 34 μL

Volume of sample B needed = 10,000 cells $\times$ 0.5 $\times$ 1.36 / 400cells/ μL = 17 μL

- 5 Calculate the sum of all of the sample volumes, V_n , to be used in the cell suspension. Using the example in **step 4**:

$$V_n = 34 \mu\text{L} + 17 \mu\text{L} = 51 \mu\text{L}$$

- 6 Calculate the volume of cold Sample Buffer, B , that is needed to bring the final volume of cell suspension to $650 \mu\text{L}$. Using the example in **step 5**:

$$B = 650 \mu\text{L} - 51 \mu\text{L} = 599 \mu\text{L}$$



For low-abundance samples, the final cell suspension can be prepared in $610 \mu\text{L}$ of cold Sample Buffer.

- 7 According to the calculations in steps 3–6, prepare the cell suspension in cold Sample Buffer in a new 1.5-mL LoBind tube.

Ensure the stock solution is well resuspended by gentle pipet-mixing before pooling.

- 8 If the samples were not filtered before counting the cells, filter through a Falcon® tube with a Cell Strainer Cap.

Loading cells in the cartridge

1 Load the cartridge with materials listed using the P1200M pipette:

Material to load	Volume (μL)	Pipette mode
Air	700	Prime/Treat
- Set P1200M pipette to Cell Load mode. - Pipet-mix the cell suspension with a manual P1000 pipette.		
Cell suspension	575	Cell Load^a
a. Press button once to aspirate 40 μL air, and then immerse tip in cell suspension. Press button again to aspirate 575 μL of cold cell suspension. Dispense 615 μL of air and cell suspension.		



Air bubbles that might appear at the inlet or outlet of the cartridge do not affect cartridge performance.

2 If necessary, wipe condensation from the top cartridge surface for optimal scanning.

3 Incubate at room temperature (15–25 °C) for 15 minutes.

During the 15-minute incubation, prepare the cell capture beads. See [Preparing BD Rhapsody™ Enhanced Cell Capture Beads](#) in the following section.



Optional: If using AbSeq, Sample Tag, or ATAC-Seq workflow, it is recommended to do two washes after the 15-minute incubation, according to the following table.

Material to load	Volume (μL)	Pipette mode
Air	700	Prime/Treat
Cold Sample Buffer	700	Prime/Treat
Air	700	Prime/Treat
Cold Sample Buffer	700	Prime/Treat

Preparing BD Rhapsody™ Enhanced Cell Capture Beads

Keep cell capture beads on ice before use.



For maximum recovery, do not vortex samples containing the beads. Gently mix suspensions with the beads by pipette only. Use low-retention pipette tips and LoBind tubes. Keep the beads cold, and pipet-mix only.

- 1 Place the beads tube on the magnet for 1 minute, and remove the storage buffer.
- 2 Remove the tube from the magnet, and pipet 750 µL of cold Sample Buffer into the tube.
- 3 Pipet-mix, and place on ice.

Loading and washing BD Rhapsody™ Enhanced Cell Capture Beads

- 1 Set the P1200M pipette to **Prime/Treat** mode.
- 2 Load the cartridge with the materials listed using the P1200M pipette:

Material to load	Volume (μL)	Pipette mode
Air	700	Prime/Treat
• Set P1200M pipette to Bead Load mode. • Use a manual P1000 to gently pipet-mix the beads in cold Sample Buffer. Immediately load.		
BD Rhapsody™ Enhanced Cell Capture Beads	630	Bead Load

- 3 Incubate the cartridge at room temperature (15–25 °C) for 3 minutes.
- 4 Place the cartridge on the plate shaker plate adapter.
- 5 Shake the cartridge at room temperature (15–25 °C) for 15 seconds.

The following is a list of suggested shakers/mixers with their corresponding settings:

- Eppendorf ThermoMixer® C: 1,000 rpm
- Eppendorf MixMate®: 1,000 rpm
- MicroPlate Genie®: 1,600 rpm. Set external timer to 15 seconds.

- 6 Blot the outlet drip with a lint-free wipe.
- 7 Return the cartridge to the Express instrument, and wait 30 seconds.
- 8 Set the P1200M pipette to **Wash** mode.
- 9 Load the cartridge with the following materials using the P1200M pipette:

Material to load	Volume (μL)	Pipette mode ^a
Air	700	Wash
Cold Sample Buffer	700	Wash
Air	700	Wash
Cold Sample Buffer	700	Wash
^a Press button once to aspirate 720 μL air or reagent. Insert the tip into the cartridge, and press button once to dispense 700 μL air or liquid. Remove pipette tip, and press button once to dispense remaining 20 μL of air or liquid.		

Lysing cells



Avoid bubbles when pipeting lysis buffer.

- 1 Add 75.0 µL of 1 M DTT to one 15-mL Lysis Buffer bottle. Check Add DTT box.



Use the Lysis Buffer with DTT ≤24 hours, and then discard.

- 2 Briefly vortex the lysis mix, and place on ice.
- 3 Move the left slider to LYSIS on the Express instrument.
- 4 Set the P1200M pipette to **Lysis** mode.
- 5 Load the cartridge with the following material using the P1200M pipette:

Material to load	Volume (µL)	Pipette mode
Lysis Buffer with DTT	550	Lysis

- 6 Incubate at room temperature (15–25 °C) for 10 minutes.



Maintain the recommended lysis time for best performance.

Retrieving BD Rhapsody™ Enhanced Cell Capture Beads

- 1 Place the 5-mL LoBind tube in the Express instrument drawer.
- 2 Ensure that the P5000M pipette is set to **Retrieval** mode.
- 3 Move the front slider to **BEADS** on the Express instrument.
- 4 Move the left slider to **RETRIEVAL**.
- 5 Leave the Retrieval magnet in the down position for 30 seconds.
- 6 Aspirate 5,000 µL of Lysis Buffer with DTT using the P5000M pipette.
- 7 Press down on the P5000M pipette to seal against the gasket.
- 8 Move the left slider to the middle position (0), and *immediately* load 4,950 µL of Lysis Buffer with DTT.
- 9 Remove the pipette from the gasket, and purge the tip.
- 10 Move the front slider to **OPEN**, and place the 5-mL LoBind tube on the large magnet with the 15-mL tube adapter for 1 minute.
- 11 Immediately proceed to [Washing BD Rhapsody™ Enhanced Cell Capture Beads](#) in the following section.
- 12 Appropriately dispose of the cartridge, Waste Collection Container, and Lysis Buffer with DTT.



All surfaces that come in contact with biological specimens can transmit potentially fatal disease. Use universal precautions when cleaning surfaces. Wear suitable protective clothing, eyewear, and gloves.

- 13 Clean the Express instrument with 10% bleach or 70% ethyl alcohol.

Washing BD Rhapsody™ Enhanced Cell Capture Beads

- 1 After the 1-minute incubation, leaving the 5-mL tube containing the retrieved beads on the large magnet, remove all but ~1 mL of supernatant without disturbing the beads.
- 2 Remove the tube from the magnet. Gently pipet-mix the beads and transfer them to a new 1.5-mL LoBind tube.
- 3 If there are still beads left in the 5-mL tube, add 0.5 mL of Lysis Buffer with DTT, rinse the 5-mL tube, and transfer to the 1.5-mL LoBind tube from the previous step.
- 4 Place the tube on the 1.5-mL magnet for ≤ 2 minutes. Remove the supernatant.



Avoid leaving Lysis Buffer or bubbles in the tube. Lysis Buffer might cause the reverse transcription reaction to fail.

- 5 Remove the tube from the magnet, and pipet 1.0 mL of cold Bead Wash Buffer into the tube. Pipet-mix.
- 6 Place the tube on the 1.5-mL magnet for ≤ 2 minutes. Remove the supernatant.
- 7 Remove the tube from the magnet, and pipet 1.0 mL of cold Bead Wash Buffer into the tube. Pipet-mix, and place on ice.



Start reverse transcription ≤ 30 minutes after washing the retrieved beads with the Bead Wash Buffer.



If performing the TCR and/or BCR assay, stop here and proceed with respective protocol to continue cDNA synthesis.

Performing reverse transcription

- 1 Ensure that the SmartBlock™ Thermoblock for ThermoMixer® C is at 37 °C.
- 2 In the pre-amplification workspace, pipet the following reagents into a new 1.5-mL LoBind tube on ice:

cDNA mix

Cap	Component	For 1 library (μL)	For 1 library + 20% overage (μL)
	RT Buffer	40.0	48.0
	dNTP	20.0	24.0
	RT 0.1 M DTT	10.0	12.0
	Bead RT/PCR Enhancer	12.0	14.4
	RNase Inhibitor	10.0	12.0
	Reverse Transcriptase	10.0	12.0
	Nuclease-Free Water	98.0	117.6
	Total	200.0	240.0

- 3 Gently vortex mix, briefly centrifuge, and place back on ice.
- 4 Place the tube of washed BD Rhapsody™ Enhanced Cell Capture Beads on the 1.5-mL tube magnet for ≤2 minutes. Remove the supernatant.
- 5 Remove the tube from the magnet and pipet 200 μL of cDNA mix into the beads. Pipet-mix.



Keep the prepared cDNA mix with beads on ice until the suspension is transferred in the next step.

- 6 Transfer the bead suspension to a new 1.5-mL LoBind tube.
- 7 Incubate the bead suspension on the thermomixer C at 1,200 rpm and 37 °C for 20 minutes.



Shaking is critical for this incubation.

- 8 Place the tube on ice.

Treating the sample with Exonuclease I

- 1 Set one thermomixer to 37 °C and a second thermomixer to 80 °C.



Exonuclease I inactivation temperatures above 80 °C can result in the loss of AbSeq molecules, thus a heat block should not be used for this step. If only one thermomixer is available, allow it to equilibrate to 80 °C before starting the inactivation incubation.

- 2 In the pre-amplification workspace, pipet the following reagents into a new 1.5-mL LoBind tube on ice:

Exonuclease I mix

Cap	Component	For 1 library (μL)	For 1 library + 20% overage (μL)
	10X Exonuclease I Buffer	20.0	24.0
	Exonuclease I	10.0	12.0
	Nuclease-Free Water	170.0	204.0
	Total	200.0	240.0

- 3 Gently vortex mix, briefly centrifuge, and place back on ice.
- 4 Place the tube of beads with cDNA mix on the 1.5-mL tube magnet for ≤2 minutes. Remove the supernatant.
- 5 Remove the tube from the magnet, and pipet 200 μL of Exonuclease I mix into the tube. Pipet-mix.
- 6 Incubate the bead suspension on the thermomixer at 1,200 rpm and 37 °C for 30 minutes.



If only one thermomixer is available, allow it to equilibrate to 80 °C before starting the inactivation incubation. Place the samples on ice until that temperature is reached.

- 7 Incubate the bead suspension on the thermomixer (no shaking) at 80 °C for 20 minutes.



Do not exceed this inactivation temperature and incubation time.

- 8 Place the tube on ice for ~1 minute.
- 9 Place the tube on the magnet for ≤1 minute until clear. Remove the supernatant.
- 10 Remove the tube from the magnet, and pipet 200 μL of cold Bead Resuspension Buffer into the tube. Pipet-mix.



Exonuclease I-treated beads can be stored at 2–8 °C for up to 1 year.

- 11 Proceed to library preparation.

Troubleshooting

For additional troubleshooting, see the troubleshooting section in the instrument user guide.

For technical support, contact your local Field Application Specialist (FAS) or scomix@bd.com.

Observation	Possible causes	Recommended solutions
No pellet after centrifuging cells or very few cells.	Rare or dilute sample.	After each centrifugation step, leave 50 µL of supernatant.

Becton, Dickinson and Company

BD Biosciences

155 North McCarthy Boulevard

Milpitas, California 95035 USA

bdbiosciences.com

scomix@bd.com