

BD Rhapsody™ Targeted mRNA and AbSeq Training Kit—4 Pack

BD Biosciences Kit Catalog No.: 633772

Purpose

The purpose of this kit is to provide the necessary reagents to support training on the BD Rhapsody™ Single-Cell Analysis system and BD Rhapsody™ Express Single-Cell Analysis system.

Contents

Box	Cat. No.	Component	Cap color	Quantity	Volume per unit	Storage
BD Rhapsody™ Cartridge Reagent Kit (Cat. No. 633731)	650000060	Cartridge Wash Buffer 1	Neutral	1 bottle	7 mL	2°C to 8°C
	650000061	Cartridge Wash Buffer 2	Neutral	1 bottle	4 mL	
	650000062	Sample Buffer	Neutral	1 bottle	28 mL	
	650000064	Lysis Buffer	Neutral	4 bottles	15 mL	
	650000065	Bead Wash Buffer	Neutral	1 bottle	10 mL	
	650000063	1 M DTT	White	1 vial	400 µL	
	650000089	Cell Capture Beads	Brown	4 vials	2 mL	
	650000090	Waste Collection Container	Neutral	4 each	—	

Box	Cat. No.	Component	Quantity	Storage
BD Rhapsody™ Cartridge Kit (Cat. No.633733)	400000847	BD Rhapsody™ Cartridge	4 each	15°C to 25°C

Box	Cat. No.	Component	Cap color	Quantity	Volume per unit	Storage
BD Rhapsody™ cDNA Kit (Cat. No. 633773)	650000076	Nuclease-Free Water	Neutral	2 vials	1 mL	-25°C to -15°C
	650000067	RT Buffer	Orange	1 vial	200 µL	
	650000068	RT 0.1M DTT	Orange	1 vial	50 µL	
	650000069	Reverse Transcriptase	Orange	1 vial	50 µL	
	650000077	dNTP	Orange	1 vial	100 µL	
	650000078	RNase Inhibitor	Orange	1 vial	50 µL	
	91-1082	Bead RT/PCR Enhancer	Black	1 vial	70 µL	
	650000071	10X Exonuclease I Buffer	Yellow	1 vial	100 µL	
	650000072	Exonuclease I	Yellow	1 vial	50 µL	
	650000066	Bead Resuspension Buffer	Black	1 vial	1 mL	

Box	Cat. No.	Component	Cap color	Quantity	Volume per unit	Storage
BD Rhapsody™ Targeted mRNA and AbSeq Amplification Kit (Cat. No. 633774)	650000076	Nuclease-Free Water	Neutral	1 vial	1 mL	-25°C to -15°C
	91-1082	Bead RT/PCR Enhancer	Black	1 vial	70 µL	
	91-1083	PCR MasterMix	White	1 vial	1.2 mL	
	91-1084	Elution Buffer	Pink	1 vial	1.8 mL	
	650000074	Universal Oligo	White	1 vial	130 µL	
	91-1085	Library Forward Primer	Red	1 vial	40 µL	
	650000080	Library Reverse Primer 1	Red	1 vial	20 µL	
	650000091	Library Reverse Primer 2	Red	1 vial	20 µL	
	650000092	Library Reverse Primer 3	Red	1 vial	20 µL	
	650000093	Library Reverse Primer 4	Red	1 vial	20 µL	
	650000066	Bead Resuspension Buffer	Black	1 vial	1 mL	
	91-1088	Sample Tag PCR1 Primer	Purple	1 vial	20 µL	
	91-1089	Sample Tag PCR2 Primer	Purple	1 vial	20 µL	
	91-1086	BD™ AbSeq Primer	Green	1 vial	70 µL	

Box	Cat. No.	Component	Cap color	Quantity	Volume per unit (µL)	Storage
BD Rhapsody™ Immune Response Panel Hs (Cat. No. 633750)	650000083	PCR1 Primers-Immune Response Panel Hs (tube label: PCR1-Immune Res. Hs)	Blue	1 vial	210	-25°C to -15°C
	650000084	PCR2 Primers-Immune Response Panel Hs (tube label: PCR2-Immune Res. Hs)	Blue	1 vial	50	

Box	Cat. No.	Component	Cap color	Quantity	Volume per unit (mL)	Storage
BD Rhapsody™ Training Cells (Cat. No. 633741)	650000081	BD Rhapsody™ Training Cells—Jurkat	Yellow	2 vials	1.0	Liquid nitrogen
	650000082	BD Rhapsody™ Training Cells—Ramos	Blue	2 vials	1.0	

Handling

Store the reagents at their specified storage temperatures. Before using reagents, see the *BD Rhapsody™ Single-Cell Analysis System Instrument User Guide* (Doc ID: 214062) or the *BD Rhapsody™ Express Single-Cell Analysis System Instrument User Guide* (Doc ID: 214063).

Safety

Use standard laboratory safety protocols. Read and understand the safety data sheets (SDSs) before handling chemicals. To obtain SDSs for chemicals ordered from BD Biosciences, go to regdocs.bd.com or contact BD Biosciences technical support at researchapplications@bd.com.

BD Biosciences

1.877.232.8995, prompt 2, 2 | researchapplications@bd.com | 2350 Qume Drive, San Jose, CA 95131 USA

Conditions: The information disclosed herein is not be construed as a recommendation to use the above product in violation of any patents. BD Biosciences will not be held responsible for patent infringement or other violations that may occur with the use of our products. Purchase does not include or carry any right to resell or transfer this product either as a stand-alone product or as a component of another product. Any use of this product other than the permitted use without the express written authorization of Becton, Dickinson and Company is strictly prohibited.

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

© 2018 BD. BD and the BD Logo are trademarks of Becton, Dickinson and Company. 07/2018

