

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

