

BD[®] Flex Single-Cell Multiplexing Kits (BD[®] Flex SMK)

A flexible sample multiplexing tool for every single-cell study on the BD Rhapsody[™] Single-Cell Analysis System



BD[®] Flex SMK



Higher sample throughput

With 24 sample tags, you can:

- Plex up to 24 samples in a single BD Rhapsody[™] Assay. Reduce sample batch effect
- Capture more than 40,000 cells per lane with <10% multiplet rate



Broader application

BD[®] Flex SMK is a species and cell type agnostic single-cell sample multiplexing tool that enables flexibility in choosing primary antibodies to tag your samples.

- More species beyond human and mouse: now you can multiplex single-cell samples from rat, chicken, pig or zebrafish in a BD Rhapsody[™] System experiment
- Labeling specific cell populations: human CD45+ cells, neurons, epithelial cells, etc.
- Multiplexing antigen-specific T-cells: anti-PE sample tags provide a seamless sample multiplexing tool for antigen-specific T-cells labeled with PE tetramers



Saving on experimental costs

- Lower sample tag cost per test or per sample with BD[®] Flex SMK
- Reduce single-cell capture and library preparation reagent cost by combining more samples to each lane. Now with 24 sample tags, your experimental expense can be lowered by up to 92% (vs no sample multiplexing)



Compatibility with all BD Rhapsody[™] Assays

BD[®] Flex SMK works seamlessly with transcriptome, protein and immune repertoire profiling assays on the BD Rhapsody[™] Single-Cell Analysis System enabling:

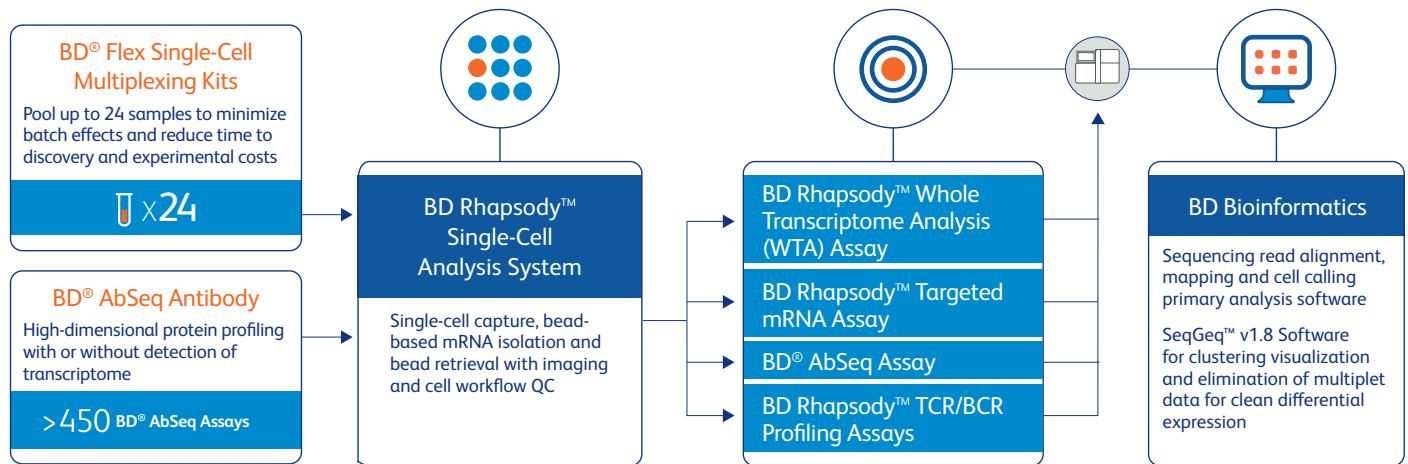
- Gene expression analysis with whole transcriptome (WTA) and targeted mRNA assays
- Cell-surface protein expression investigation with or without transcriptome detection using BD[®] AbSeq Assays. BD offers more than 450 individual AbSeq antibody-oligonucleotides and 30-plex BD[®] AbSeq Immune Discovery Panel. Custom BD[®] AbSeq Antibody-Oligonucleotides are also available upon request
- Immune repertoire discovery with full-length B-cell and T-cell receptor profiling assays or V(D)J CDR3 protocols



Reliable sample multiplexing performance

- Sample tags optimized to test sizes, validated for 50K–1M cell staining per test
- Comparable sensitivity and selectivity as one-step antibody-oligo labeling
- Multiplet identification for cleaner single-cell sequencing analysis





Ordering information

Description	Cat. No.
BD® Flex Single-Cell Multiplexing Kit A	633849
BD® Flex Single-Cell Multiplexing Kit B	633850
BD® Flex Single-Cell Multiplexing Kit C	633851
BD® Flex Single-Cell Multiplexing Kit D	633852

To request a quote or place an order, visit bdbiosciences.com or contact your local BD sales representative.

Suggested companion products

Description	Cat. No.
BD® AbSeq Immune Discovery Panel	Contact for more info
BD® AbSeq Single-Vial Reagents	Contact for more info
BD Pharmingen® Human BD Fc Block® Buffer Solution	564219
BD Pharmingen® Purified Rat Anti-Mouse CD16/CD32 (Mouse BD Fc Block®)	553141
BD Pharmingen® Purified Mouse Anti-Rat CD32 (Rat BD Fc Block®)	550270
BD Rhapsody™ Whole Transcriptome Analysis (WTA) Amplification Kit	633801
BD Rhapsody™ Targeted mRNA and AbSeq Amplification Kit	633774
BD Rhapsody™ cDNA Kit	633773
BD Rhapsody™ Enhanced Cartridge Reagent Kit	664887
BD Rhapsody™ 8-Lane Cartridge	666262
BD Rhapsody™ HT Xpress Package	666625
BD Rhapsody™ Single-Cell Analysis System	633701

To request a quote or place an order, email scomix@bdscomix.com or contact your local BD sales representative.

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

BD Life Sciences, San Jose, CA 95131, USA

bdbiosciences.com

