

Data Sheet

BCMA, Fc-Fusion, Avi-Tag, PE-Labeled

Human, Recombinant, C-terminal Fc-Avi-Tag,
PE-Labeled

Catalog #: 100733

Lot#: 200408

Conc.: 1.97 mg/ml

Formulated in: 8 mM Phosphate, pH 7.4,
110 mM NaCl, 2.2 mM KCl, and 20%
glycerol.

Stability: At least 6 months at -80°C . *Avoid
freeze/thaw cycles.*

References:

1. Schiemann, B., et al., Science. 2001;
293(5537): 2111-2114.
2. Gross, J., et al., Nature. 2000;
404:995-999.

Description: Human BCMA, known as TNFRSF17, CD269, and TNFRSF13A, GenBank Accession No. NM_001192, a.a. 1-54, with C-terminal Avi-tag fused to the Fc portion of Human IgG1, and labeled with phycoerythrin (PE), expressed in a HEK293 cell expression system. MW = 34 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Applications: Useful for labeling cells expressing Anti-BCMA antibodies for flow cytometry and immunofluorescence microscopy.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:

8 μg BCMA-RPE

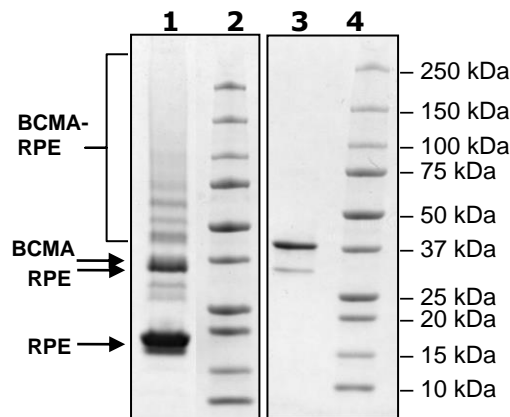
Lane 3:

4 μg BCMA

Lane 2 and 4:
Protein Marker

MW: BCMA 34 kDa

Purity: $\geq 90\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

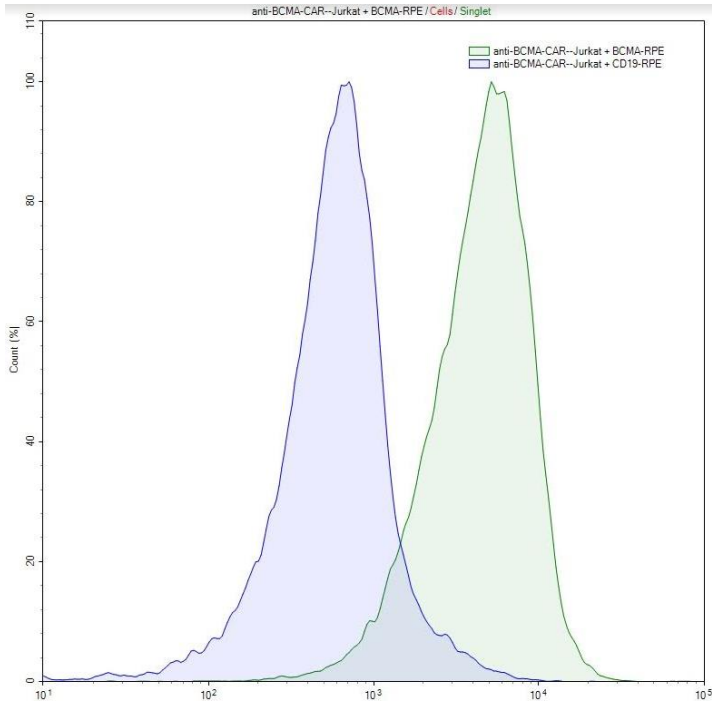
To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com



BCMA, Fc-Fusion, Avi-Tag, PE-Labeled (BPS Bioscience #100733) was tested using FACS with Anti-BCMA CAR /NFAT (Luciferase) Reporter Jurkat Cell Line (BPS Bioscience #79694) and Anti-CD19 CAR / NFAT (Luciferase) Reporter Jurkat Cell Line (CD19 SCFV-CD28-4-1BB-CD3 ζ) (BPS Bioscience #79853).

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com