

Data Sheet

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

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

Catalog #: 79467

Lot#: 190124-1

Conc.: 0.50 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, BCM, BCMA, 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 expressed in a HEK293 cell expression system. This protein runs at a higher MW by SDS-PAGE due to glycosylation. This protein is enzymatically biotinylated using Avi-Tag™ technology. MW = 34 kDa.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:

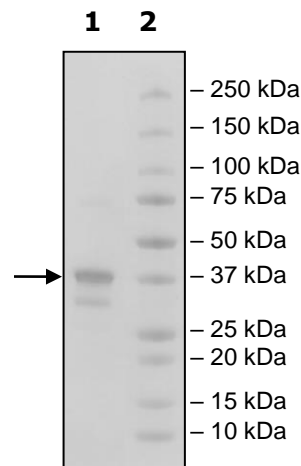
4µg BCMA

Lane 2:

Protein Marker

MW: 34 kDa + glycans

Purity: $\geq 73\%$



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