



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

Data Sheet

Anti-BCMA-Anti-CD19-Anti-CD3-His Trispecific Antibody

Catalog #:100761

Lot #: 200513 Conc.: 0.24 mg/ml

Description: The Anti-BCMA-Anti-CD19-Anti-CD3-His Trispecific Antibody is engineered to bind to three different targets simultaneously: 1) BCMA, a B cell antigen that is highly expressed by mature B cells and malignant myeloma cells, 2) CD19, a B-cell marker that is a target for many leukemias and lymphomas, and 3) CD3, a T cell receptor activation signal that leads to cytokine secretion. The multi-functionality of this trispecific antibody allows it to bind to BCMA and CD19 on the tumor cell and CD3 on T cells simultaneously, thus bringing T lymphocytes closer to the cancer cells. The binding event targets the tumor while providing co-stimulatory signals that promote T cell expansion and cytotoxicity against BCMA+ and CD19+ cancer cells.

Background: B-cell maturation antigen (BCMA), also known as tumor necrosis factor receptor superfamily member 17 (TNFRSF17), is a protein encoded by the TNFRSF17 gene. TNFRSF17 is a cell surface receptor of the TNF receptor superfamily that recognizes B-cell activating factor (BAFF). BCMA is preferentially expressed in mature B lymphocytes and also on Multiple Myeloma (MM) cells. Upregulation of BCMA also correlates with disease burden and prognosis in multiple myeloma.

B-lymphocyte antigen CD19 (Cluster of Differentiation 19), also known as B-Lymphocyte Surface Antigen B4 and CVID3, is a transmembrane protein expressed in follicular dendritic cells and all B lineage cells except plasma cells. CD19 plays two major roles in human B cells. It acts as an adaptor protein to recruit cytoplasmic signaling proteins to the membrane and it works within the CD19/CD21 complex to decrease the threshold for B cell receptor signaling pathways. Due to its presence on all B cells, it is a biomarker for B lymphocyte development and lymphoma diagnosis and can be used as a target for leukemia immunotherapies.

Application: This product is for research use only. It is not suitable for human diagnostic or therapeutic use. The anti-BCMA-anti CD19-anti-CD3 can be used for studying BCMA+ cancer cell-mediated T cell activation, using either primary T cells or reporter cell lines such as NFAT-luc-Jurkat cells (BPS Bioscience #60621).

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

Host Species: Human

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

Purification: His-tag affinity purification from HEK293 cells

Stability: Stable for at least 12 months at -80°C. Avoid freeze/thaw cycles.

Fig. 1: Protein expression of Anti-BCMA-anti-CD19-anti-CD3-His-Tag antibody

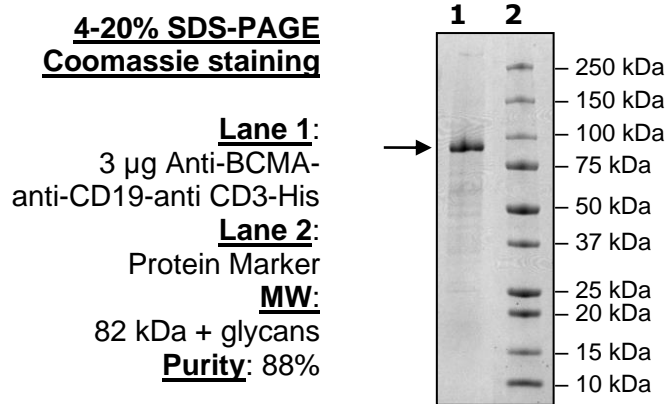
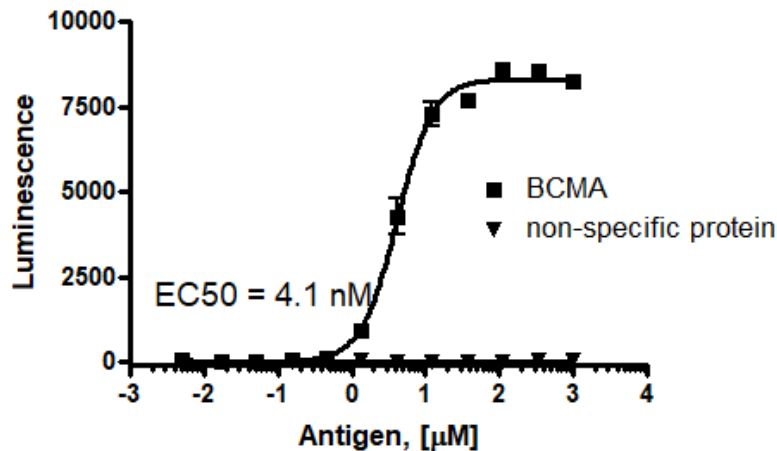


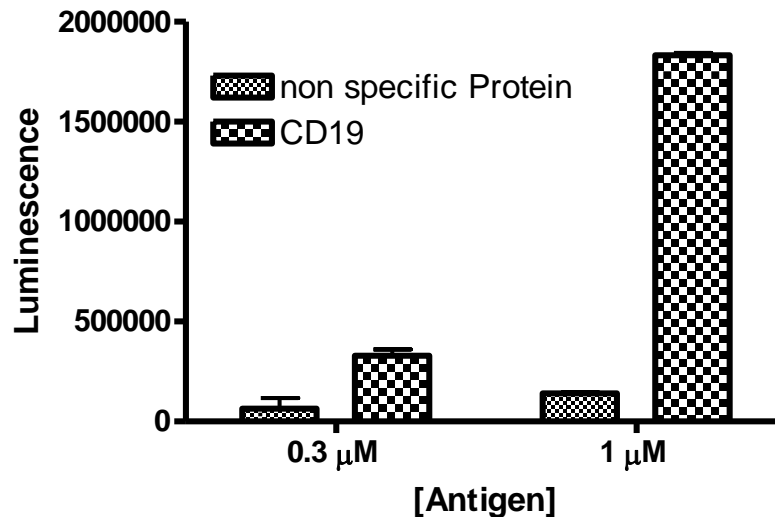
Fig. 2: Binding Specificity of anti-BCMA-anti-CD19-anti-CD3-His antibody to BCMA. Titration of BCMA Biotin (BPS Bioscience #79467-1) or control non-specific biotinylated protein binding to anti-BCMA-anti-CD19-anti-CD3-His coated on an ELISA plate, followed by streptavidin-HRP and chemiluminescent detection.



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

Fig. 3: Binding Specificity of anti-BCMA-anti-CD19-anti-CD3-His antibody to CD19. Binding of CD19-Biotin (BPS Bioscience #79475) or control non-specific biotinylated protein to anti-BCMA-anti-CD19-anti-CD3-His coated on an ELISA plate, followed by streptavidin-HRP and chemiluminescent detection.



Experimental design and assay protocol for measuring anti-BCMA-anti-CD19-anti-CD3 antibody functional activity using NFAT-luc reporter Jurkat cell line:

Jurkat effector cells with endogenous TCR/CD3 and transfected reporter NFAT-luc (BPS Bioscience #60621) are incubated with increasing concentrations of anti-BCMA-anti-CD19-anti-CD3 antibody in the presence of BCMA-CHO cells (BPS Bioscience #79500-H), CD19-CHO cells (BPS Bioscience #79561-H), or parental CHO cells (ATCC #CCL-61™).

1. Seed CHO, BCMA-CHO or CD19-CHO cells at 30,000 cells/well in a 96-well clear bottom white plate and allow a few hours for the cells to attach.
2. Add 30,000 NFAT-luc Jurkat cells/well.
3. Add the antibody at a recommended dilution range of 100 fM-100 nM. The tri-specific antibody simultaneously binds to TCR/CD3 on Jurkat/NFAT reporter cells and tumor antigen BCMA on BCMA-CHO cells or CD19 on CD19-CHO cells. The total volume of CHO cells, NFAT-luc Jurkat cells, and the antibody was adjusted to 125ul per well.

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

- After 16 hours the luciferase activity is measured by ONE-Step™ luciferase assay (BPS Bioscience #60690) per recommended protocol. The bispecific antibody interaction stimulates NFAT-luciferase activity.

Fig. 4: Activation of NFAT Reporter Jurkat cells by anti-BCMA-anti-CD19-anti-CD3 antibody in the presence of BCMA-CHO cells or CHO cells. EC50=62.9 pM.

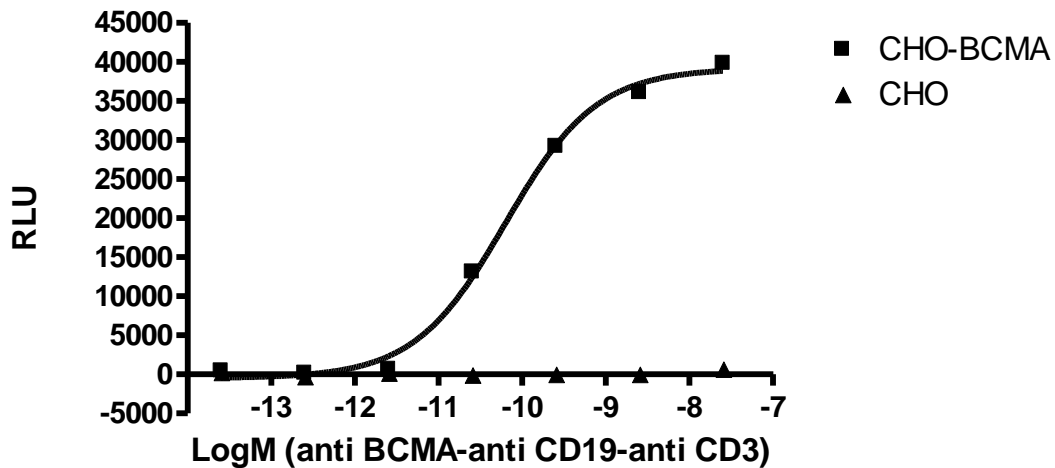
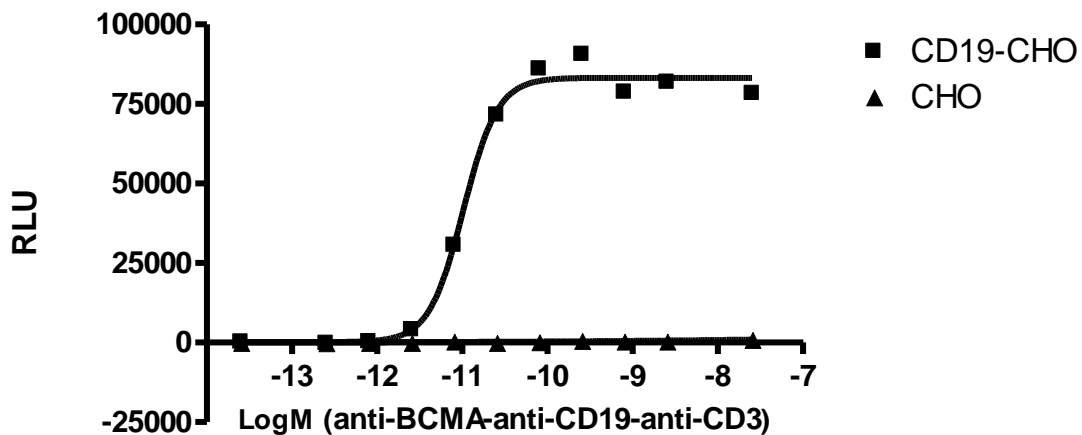


Fig. 5: Activation of NFAT Reporter Jurkat cells by anti-BCMA-anti-CD19-anti-CD3 antibody in the presence of CD19-CHO cells or CHO cells. EC50=10.3 pM.



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

Related Products:

<u>Product</u>	<u>Cat. #</u>	<u>Size</u>
NFAT Reporter (Luc) – Jurkat Cell Line	60621	2 vials
BCMA CHO Cell Line (High Expression)	79500-H	2 vials
CD19 CHO Cell Line (High Expression)	79561-H	2 vials
ONE-Step™ Luciferase Assay System	60690-1	10 ml
BCMA, Fc-fusion (IgG1), Avi-Tag, Biotin-Labeled	79467-1	25 µg
CD19, Fc-fusion (IgG1), Avi-Tag, Biotin-Labeled	79475-1	25µg
Anti-BCMA Antibody (Single-Chain Variable Fragment)	100173	50 µg
Anti-CD19-Anti-CD3 Bispecific Antibody	100441	50 µg
Anti-BCMA-Anti-CD3 Bispecific Antibody	100689	50 µg

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