

Data Sheet

Anti-BCMA Antibody (Single-Chain Variable Fragment), His-Tag

Recombinant, Human, Monoclonal,
C-terminal His-Tag

Catalog #: 100173

Lot #: 180815-2

Conc.: 1.96 mg/ml

Host Species | Isotype:

Human | IgG1

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

Stability: Stable for at least 12 months at
-80°C. *Avoid freeze/thaw cycles. Store
diluted antibody with carrier protein (BSA 0.1
– 0.5%) at 4°C, stable for 2 weeks.*

References:

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

Description: Single-chain variable fragment
Anti-BCMA recombinant human antibody with
C-terminal His-Tag. This antibody not been
tested for cross-reactivity with other species.

Purification: Protein A affinity
chromatography from HEK293 supernatants.

Quality Assurance

Anti-BCMA: BCMA (Biotin) Binding

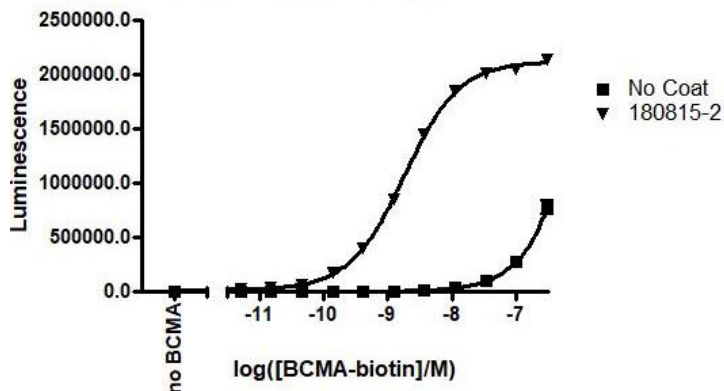


Fig.1 Antibody coat proteins were added at 100ng/well and Biotin-labeled BCMA (BPS #79467) was titrated from 0-300 nM with a 3-fold dilution series. Binding was detected with Strep-HRP.

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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com