

Data Sheet

Spike S1 Neutralizing Antibody (VHH), Fc-fusion (IgG1), Avi-Tag (SARS-CoV-1), Biotin-labeled

SARS-CoV-1, Monoclonal, C-terminal Fc-Avi-Tag
Catalog #: 100785
Lot #: 200604 **Conc.:** 1.00 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 6 months at -80°C.
 Avoid freeze/thaw cycles.

References:

1. Ho, M. *et al.*, *Antibody Therapeutics*. 2020; **3(2)**: 109-114
2. Wrapp, D. *et al.*, 2020. *Cell* **181**: 1004-1015

Description:

SARS-CoV-1 monoclonal Spike S1 VHH neutralizing antibody, a.a. 1-127(full length), with C-terminal Avi-Tag fused to the Fc portion of Human IgG1, expressed in a HEK293 cell expression system. MW = 43 kDa. This protein runs at a higher M.W. by SDS-PAGE due to glycosylation. This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%. VHHs are single-domain antibodies from camelids.

Purification:

Protein A Chromatography.

Application:

ELISA and neutralization assays.

Quality Assurance

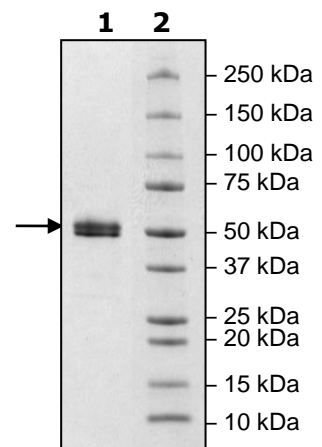
**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 4µg SARS-CoV-1

Lane 2:
 Protein Marker

MW: 43 kDa+glycans

Purity: ≥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

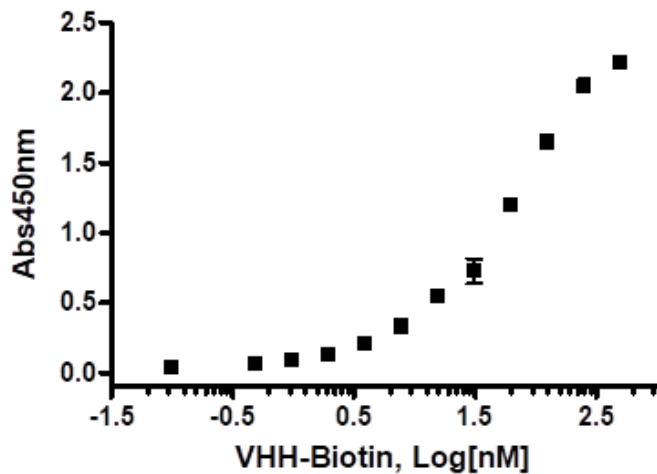


Fig.1 Spike Trimer (S1+S2), His-tag (SARS-CoV-1) (BPS Bioscience, #100789) and Spike S1 Neutralizing Antibody (VHH), Fc-fusion (IgG1), Avi-Tag (SARS-CoV-1), Biotin-labeled binding was measured by ELISA using Streptavidin-HRP and Colorimetric HRP substrate.

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