

Data Sheet

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

SARS-CoV-1, Monoclonal, C-terminal Fc-Avi-Tag
Catalog #: 100784
Lot #: 200603 **Conc.:** 1.35 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. VHHs are single-domain antibodies from camelids.

Purification:

Protein A Chromatography.

Application:

ELISA and neutralization assays.

Quality Assurance

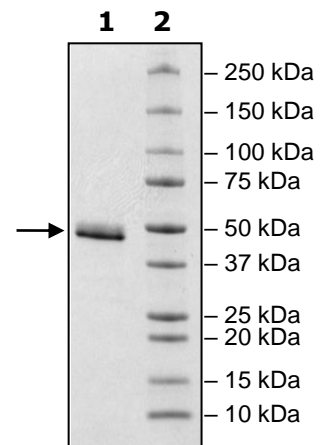
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
4µg SARS-CoV-1
(VHH)

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

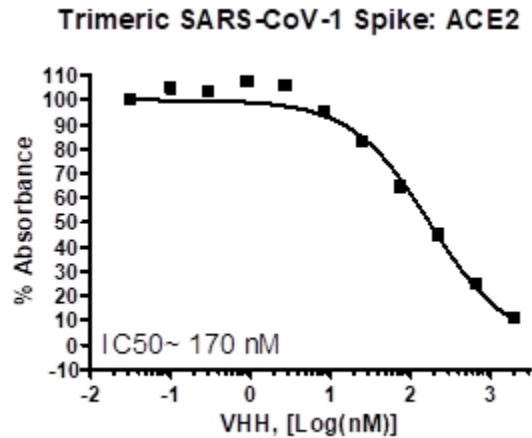


Fig.1 Inhibition of ACE2:SARS-CoV-1 Spike binding by a small anti-SARS-CoV-1 Spike antibody, VHH (BPS Bioscience, #100784). A small anti-SARS-CoV-1 Spike antibody was evaluated using the SARS-CoV-1 Spike Trimer (S1+S2):ACE2 Inhibitor Screening Colorimetric Assay Kit (BPS Bioscience, #79999). The antibody was serially diluted in 1x Immuno Buffer (BPS Bioscience, #79311) and tested following the assay kit protocol.

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