

## Data Sheet

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

SARS-CoV, 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 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

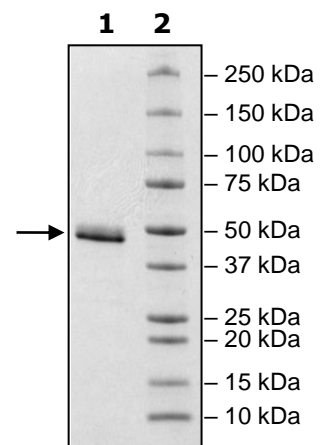
**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
4µg SARS-CoV  
(VHH)

**Lane 2:**  
Protein Marker

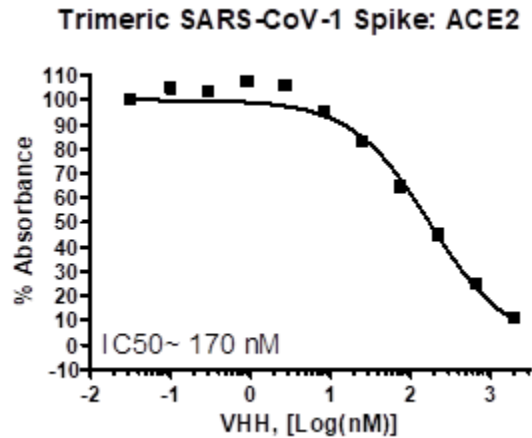
**MW:** 43 kDa+glycans

**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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)