

Data Sheet

Spike S1, Fc Fusion, Avi-tag (SARS-CoV-2)

SARS-CoV-2, Recombinant, C-terminal
 Fc-Avi-tag
Catalog #: 100678
Lot #: 200325 **Conc:** 0.57 mg/ml

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

Stability: At least 6 months at -80°C. *Avoid
 freeze/thaw cycles.*

References:

1. Hoffman, M. *et al.*, *Cell*. 2020 Apr;
181:1-10.
2. Walls, A. C. *et al.*, *Cell*. 2020 Mar;
180:1-12

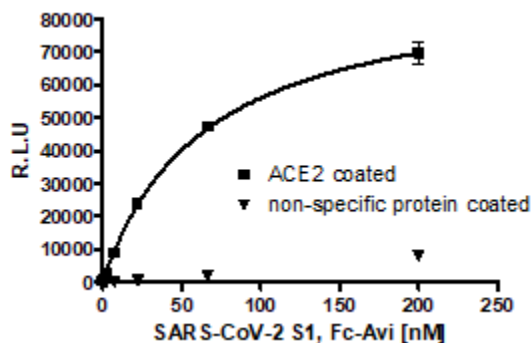
Description: Severe acute respiratory
 Coronavirus 2 Spike Glycoprotein S1 (SARS-
 CoV-2 Spike S1), GenBank Accession No.
 QHD43416.1, a.a. 13-665, fused at the C-
 terminus of the Fc portion of human IgG1, with
 a C-terminal Avi-tag™, expressed in a
 HEK293 expression system, MW=102 kDa.
 This protein runs at a higher MW by SDS-
 PAGE due to glycosylation.

Assay Conditions: SARS-CoV-2 S1 and
 ACE2 binding assay was done using BPS
 ACE2:SARS-CoV-2 Spike Inhibitor Screening
 Assay Kit (BPS Bioscience #79936). SARS-
 CoV-2 S1 protein and anti-human Fc-HRP
 were used instead of SARS-CoV-2 Spike
 protein (RBD) and anti-mouse Fc-HRP
 respectively. Non-specific His-tag protein was
 coated and used for a negative control.

Applications: Useful for the study of enzyme
 kinetics, screening inhibitors, and selectivity
 profiling.

Quality Assurance

SARS-CoV-2 S1:ACE2 Binding

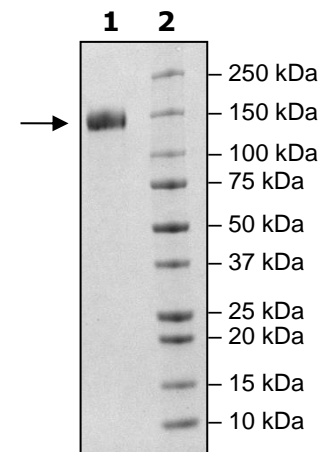


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
 6µg SARS-CoV-2 S1
Lane 2:
 Protein Marker

MW: 102 kDa

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