

Data Sheet

Spike S1, Fc fusion, Avi-tag, Biotin-Labeled (SARS-CoV-2)

SARS-CoV-2, Recombinant, C-terminal
 Fc-Avi-tag, Biotin-labeled
Catalog #: 100679
Lot #: 200326 **Conc:** 0.37 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

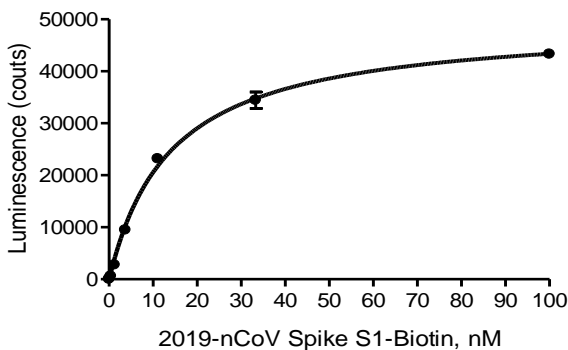
Description: Severe acute respiratory Coronavirus 2 Spike Glycoprotein S1 (SARS-CoV-2 Spike S1), GenBank Accession No. QHD43416.1, a.a. 13-665, fused at the C-terminus of the Fc portion of human IgG1, with a C-terminal Avi-tag, expressed in a HEK293 expression system and enzymatically biotinylated using Avi-tag™ technology. Biotinylation is confirmed to be ≥90%. MW=102 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Assay Conditions: Nickel plate wells were coated with 1 µg/ml ACE2, His-Tag (BPS Bioscience, #11003, 50 ng/well). Then, 2019-nCoV Spike S1, Fc-fusion (IgG1), Biotin-labeled was incubated with ACE2, His-tag for 1 hour at room temperature, 50 µl/well. Bound protein was quantified with Streptavidin-HRP (BPS Bioscience, #79742) and luminescence was measured in a BioTek Synergy 2 microplate reader.

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

Quality Assurance

ACE2:SARS-CoV-2 Spike Interaction



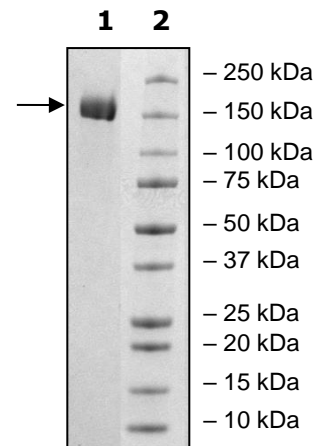
4-20% SDS-PAGE Coomassie staining

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