

## 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 #:** 100720  
**Lot #:** 200423                      **Conc:** 1.02 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*. 2002 Mar;  
**180**:1-12

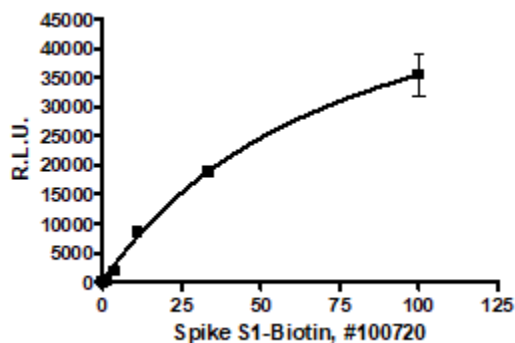
**Description:** Severe acute respiratory  
Coronavirus 2 Spike Glycoprotein S1 (SARS-CoV-  
2 Spike S1), GenBank Accession No.  
QHD43416.1, a.a. 16-685, fused at the C-terminus  
of the Fc portion of human IgG1, with a C-terminal  
Avi-tag™, expressed in a HEK293 expression  
system enzymatically biotinylated using Avi-tag™  
technology. Biotinylation is confirmed to be ≥90%.  
MW=104 kDa. This protein runs at a higher  
molecular weight due to glycosylation.

**Assay Conditions:** SARS-CoV-2 S1(BPS  
Bioscience #100720) and ACE2 binding assay  
was done using BPS ACE2:SARS-CoV-2 Spike  
Inhibitor Screening Assay Kit (BPS Bioscience  
#79936). SARS-CoV-2 S1 (BPS Bioscience  
#100720) protein and Streptavidin-HRP were  
used instead of SARS-CoV-2 Spike protein (RBD)  
and anti-mouse Fc-HRP respectively.

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

### Quality Assurance

#### ACE2:SARS-CoV-2 Inhibition

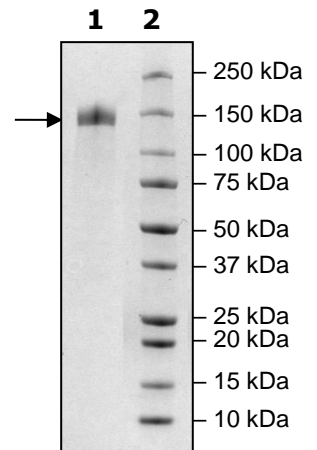


#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
4µg SARS-CoV-2 S1  
**Lane 2:**  
Protein Marker

**MW:** 104kDa

**Purity:** ≥90%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

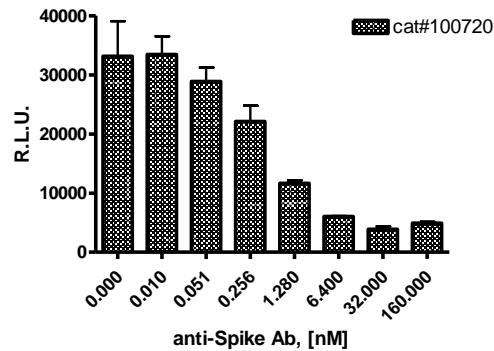
To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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)



6042 Cornerstone Court West, Suite B  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)



SARS-CoV-2 S1(BPS Bioscience #100720, 20 nM) was incubated with various concentrations of the anti-Spike Ab for 1 hour at room temperature. After then, S1/Ab mixture was added to the plate coated with ACE2 as described in the assay kit (BPS Bioscience #79936). Streptavidin-HRP were used to measure SARS-CoV-2-S1/ACE2 binding.

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)