

## Data Sheet

### Spike S1, Mouse Fc-tag (SARS-CoV-2)

SARS-CoV-2, Recombinant,  
C-terminal mFc-tag

**Catalog #:** 100684

**Lot #:** 200506

**Conc.:** 0.55 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  $-80^{\circ}\text{C}$ . *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**(2):281-292.

**Description:** Severe acute respiratory Coronavirus 2 Spike Glycoprotein S1 (SARS-CoV-2 Spike S1), also known as novel coronavirus spike S1 and nCoV spike S1, GenBank Accession No. QHD43416.1, a.a. 319-541, with a C-terminal mouse Fc-tag (mFc), expressed in a HEK293 cell expression system. MW=50 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

## Quality Assurance

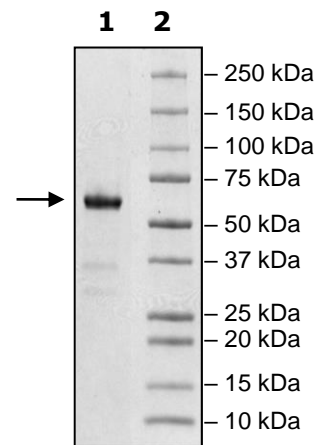
### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
4 $\mu\text{g}$  SARS-CoV-2(mFc)

**Lane 2:**  
Protein Marker

**MW:** 50 kDa

**Purity:**  $\geq 90\%$



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)