

## Data Sheet

### 2019-nCoV Spike S1, Fc-tag

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

**Catalog #:** 100688

**Lot #:** 200325

**Conc.:** 0.05 mg/ml

**Formulated in:** 50 mM Tris-HCl, pH 7.5, 90 mM glycine, 0.1 mM PMSF.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

#### References:

1. Zhou, P. *et. al*, *Nature*. **272**: 270-289, 2020.
2. Xiao, X. *et. al*, *Cell Mol Life Sci*. **61(19-20)**: 2428-2430, 2004.

**Description:** SARS-CoV-2 2019-nCoV Spike protein S1, also known as SARS-CoV s1 and coronavirus spike S1, GenBank Accession No. QHD43416.1, a.a. 16-685, with C-terminal Fc-tag, expressed in a CHO cell expression system. MW= 160 kDa.

## Quality Assurance

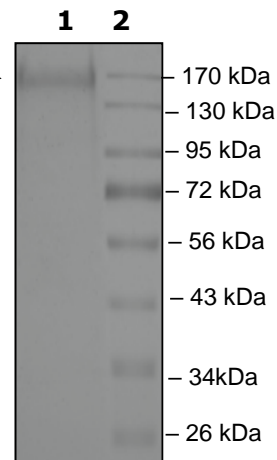
#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:** →  
2019-nCoV Spike S1

**Lane 2:**  
Protein Marker

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

**MW:** 160 kDa



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)