

Data Sheet

Spike S2, Fc-Tag **(SARS-CoV-2)**

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

Catalog #: 100895

Lot #: 200917

Conc.: 0.40 mg/ml

Description: SARS-CoV-2 Spike protein S2 subunit, also known as 2019-nCoV Spike S2, GenBank Accession No. MN908947, a.a. 686-1212, with C-terminal Fc-tag, expressed in a CHO cell expression system. MW=130 kDa.

Formulated in: 50 mM Tris-HCl, pH 7.5, 90 mM Glycine.

Stability: At least 6 months at -80°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.

Quality Assurance

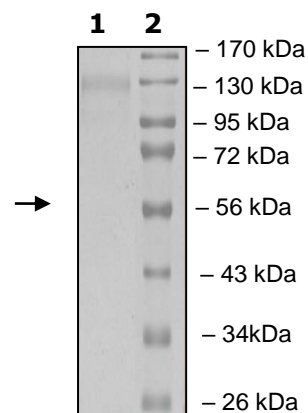
4-20% SDS-PAGE **Coomassie staining**

Lane 1:
Spike S2

Lane 2:
Protein Marker

Purity: >90%

MW: 130 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

Please visit our website at: www.bpsbioscience.com