

Data Sheet

Spike S1, Avi-His-tag (SARS-CoV-2)

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

Catalog #: 100730

Lot #: 200429

Conc: 2.13 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(2):281-292.

Description: Severe acute respiratory
Coronavirus 2 Spike Glycoprotein S1 (SARS-
CoV-2 Spike S1), GenBank Accession No.
QHD43416.1, a.a. 16-685 with a C-terminal
Avi-His-tag, expressed in a HEK293
expression system, MW=78 kDa. This protein
runs at a higher MW by SDS-PAGE due to
glycosylation.

Quality Assurance

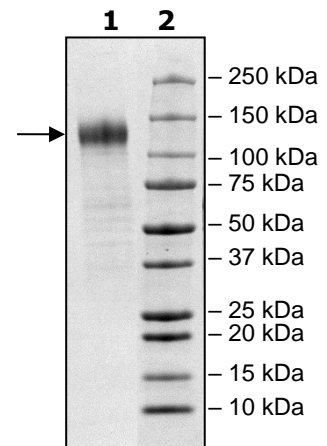
4-20% SDS-PAGE Coomassie staining

Lane 1:
4µg SARS-CoV-2 S1

Lane 2:
Protein Marker

MW: 78 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