

Data Sheet

Spike Trimer (S1+S2), His-tag (SARS-CoV)

SARS-CoV, Recombinant, C-terminal
His-tag

Catalog #: 100789

Lot #: 200610

Conc: 0.75 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.

Description: Severe acute respiratory Coronavirus SARS Coronavirus Spike trimer (S1+S2) (SARS-CoV S protein), Genbank Accession No. AAP13567, a.a. 1-1195(full length), with a C-terminal His-tag, expressed in a HEK293 expression system. MW=136 kDa. This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.

References:

1. Watanbae, Y. *et al.*, *Science*. 2020 May; DOI: 10.1126/science.abb9983
2. Jaume, M. *et al.*, *J. Virology*. 2011 Oct; **85(20)**: 10582-10597

Quality Assurance

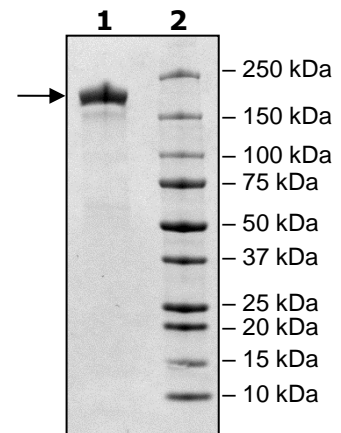
4-20% SDS-PAGE Coomassie staining

Lane 1:
4µg SARS-CoV1 S
Trimer

Lane 2:
Protein Marker

MW: 136 kDa+glycans

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