

Data Sheet

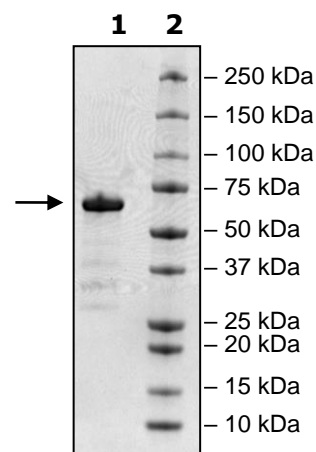
Spike S1 RBD, Fc-Fusion, Avi-Tag (SARS-CoV-2)

SARS-CoV-2, Recombinant,
C-terminal Fc-Avi-tag**Catalog #:** 100698**Lot #:** 200429**Conc.:** 1.43 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. 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 Spike protein S1 subunit, receptor binding domain (RBD), also known as SARS-CoV-2 spike RBD, novel coronavirus spike RBD and nCoV spike RBD, GenBank Accession No. MN_908947.1, a.a. 319-541, fused at the C-terminus of the Fc portion of human IgG1, with a C-terminal Avi-tag™, expressed in a HEK293 cell expression system. MW=54 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 μg SARS-CoV-2
RBD**Lane 2:**
Protein Marker**MW:** 54 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.comPlease visit our website at: www.bpsbioscience.com