

Data Sheet

Spike S1 RBD (V367F), Avi-His-tag **Biotin-labeled (SARS-CoV-2)**

SARS-CoV-2, Recombinant, C-terminal
Avi-His-tag, Biotin-labeled

Catalog #: 100770

Lot #: 200529-2

Conc.: 1.15 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 S1 receptor binding domain (RBD), also known as SARS-CoV-2 Spike 1 RBD, novel coronavirus Spike 1 RBD and nCoV Spike 1 RBD, GenBank Accession No. QHD43416.1, a.a. 319-541 with a V367F mutation and a with C-terminal Avi-His-tag, expressed in a HEK293 cell expression system and enzymatically biotinylated using Avi-tag™ technology. Biotinylation is confirmed to be $\geq 90\%$. MW=28 kDa. This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.

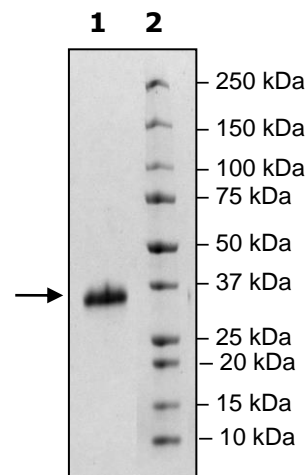
Quality Assurance

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

Lane 1:
4 μg SARS
RBD(V367F)
Lane 2:
Protein Marker

MW: 28 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**

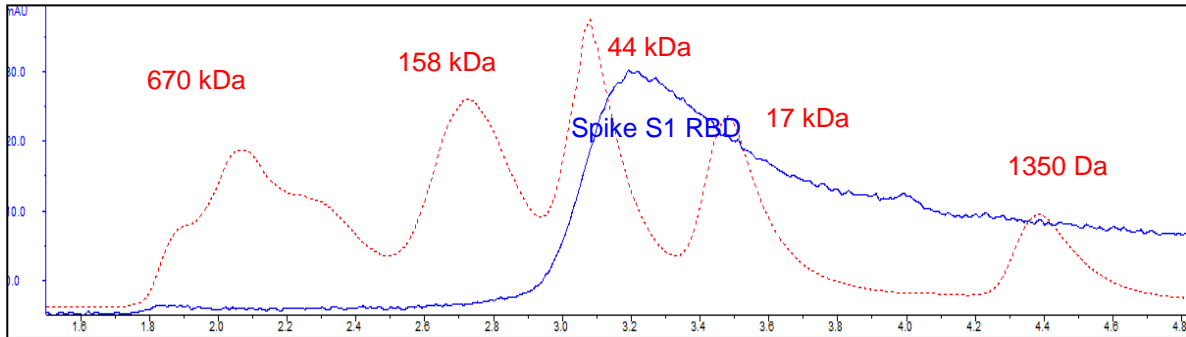
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Gel Filtration



<10% Aggregation

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