

## Data Sheet

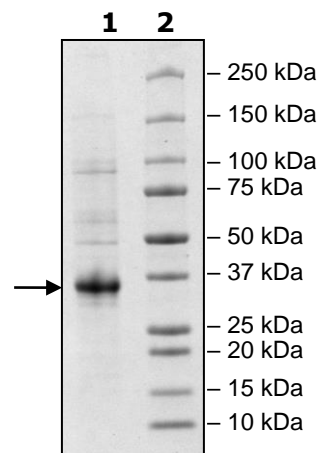
### **Spike S1 RBD, Avi-His-tag, Biotin-labeled (SARS-CoV-2)**

SARS-CoV-2, Recombinant, C-terminal  
Av-His-tag, Biotin-labeled**Catalog #:** 100697**Lot #:** 200529**Conc.:** 0.70 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  $-80^{\circ}\text{C}$ . *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 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:**  
5 $\mu\text{g}$  SARS-CoV-2  
RBD**Lane 2:**  
Protein Marker**MW:** 28 kDa**Purity:**  $\geq 79\%$ 

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](mailto:info@bpsbioscience.com)Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)