

## Data Sheet

### **Spike S1 RBD (G476S), Avi-His-tag** **(SARS-CoV-2)**

SARS-CoV-2, Recombinant, C-terminal  
Avi-His-tag**Catalog #:** 100868**Lot #:** 200902**Conc.:** 0.62 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:**

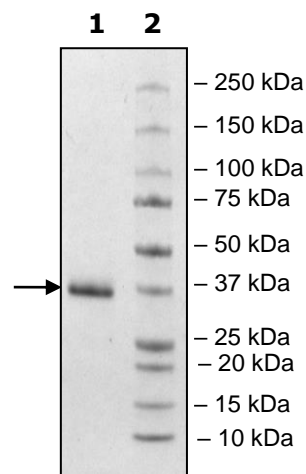
1. Zhou, P. *et al.*, *Nature*. **272**: 270-289, 2020.
2. Xiao, X. *et al.*, *Cell Mol Life Sci*. **61(19-20)**: 2428-2430, 2004.

**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 G476S mutation and with a C-terminal Avi-His-tag, expressed in a HEK293 cell expression system. 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:**3 $\mu\text{g}$  SARS  
RBD(G476S)**Lane 2:**

Protein Marker

**MW:** 28 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](mailto:info@bpsbioscience.com)Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)