

Data Sheet

2019-nCoV Spike RBD, His-tag

SARS-CoV-2, Recombinant, C-terminal His-tag
Catalog #: 100687
Lot #: 200325 **Conc.:** 0.20 mg/ml

Formulated in: 50 mM sodium phosphate, pH 7.5, 300m M NaCl, 150 mM imidazole.

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.

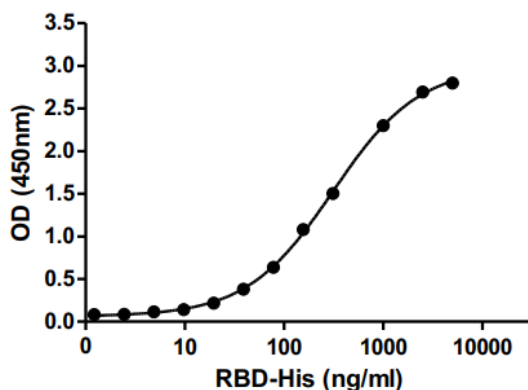
Description: SARS-CoV-2 2019-nCoV Spike protein S1 subunit, receptor binding domain (RBD), also known as SARS-CoV-2 spike RBD, novel coronavirus spike RBD and nCoV spike RBD, GenBank Accession No. QHD43416.1, a.a. 319-541, with C-terminal His-tag, expressed in a CHO cell expression system. MW= 39 kDa.

Assay Conditions: Measured by its binding ability in a functional ELISA, immobilized human ACE2, Fc-tag protein (a.a. 19-740) at 2 µg/ml (100 µl/well) can bind to 2019-nCoV spike protein RBD, His-tag.

Applications: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

ACE2:2019-nCoV Spike RBD Binding



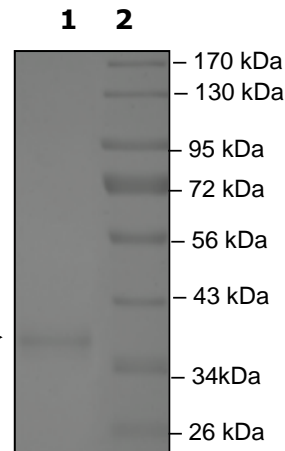
4-20% SDS-PAGE Coomassie staining

Lane 1:
2019-nCoV Spike RBD

Lane 2:
Protein Marker

Purity: ≥90% →

MW: 39 kDa



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