

Data Sheet

Spike S1 RBD-Nucleocapsid Protein Chimera (SARS-CoV-2)

SARS-CoV-2, Recombinant, Chimera

Catalog #: 100938

Lot #: 220405

Conc.: 0.20 mg/ml

Formulated in: 1x PBS, pH 7.4

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

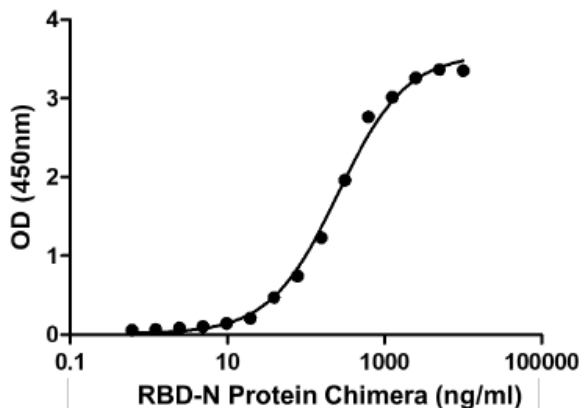
References:

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 (Spike S1 RBD), GenBank Accession No. MN908947, a.a. 319-541, fused with HSA to SARS-CoV-2 Nucleocapsid protein (N-protein), GenBank Accession No. QHD43423, a.a. 237-419, with C-terminal His-tag, Expressed in CHO cells. MW=122 kDa.

Quality Assurance

Binding of Spike S1 RBD-Nucleocapsid Protein Chimera (SARS-CoV-2) to Nucleocapsid Protein hlgG antibody (SARS-CoV-2)



Binding of Spike S1 RBD-Nucleocapsid Protein Chimera (SARS-CoV-2) to immobilized Nucleocapsid Protein hlgG antibody (SARS-CoV-2) determined by ELISA using HRP-conjugated Anti-Spike hlgG antibody (SARS-CoV-2) as a detection antibody.

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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com