

Data Sheet

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

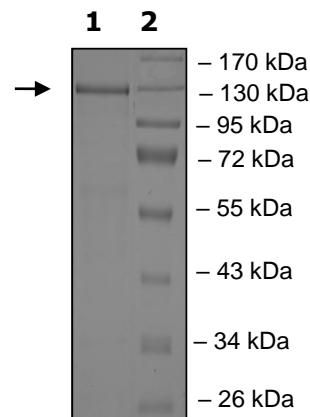
SARS-CoV-2, Recombinant, Chimera

Catalog #: 100938**Lot #:** 201027**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=130 kDa.**Applications:** Used for the detection and quantification of antibodies to SARS-CoV-2 spike protein S1 (RBD) and Nucleocapsid (N) Protein from biological samples.

Quality Assurance

**4-20% SDS-PAGE
Coomassie staining****Lane 1:
Spike S1 RBD-
Nucleocapsid
Protein Chimera****Lane 2:
Protein Marker****Purity: $\geq 70\%$** **MW: 130 kDa**

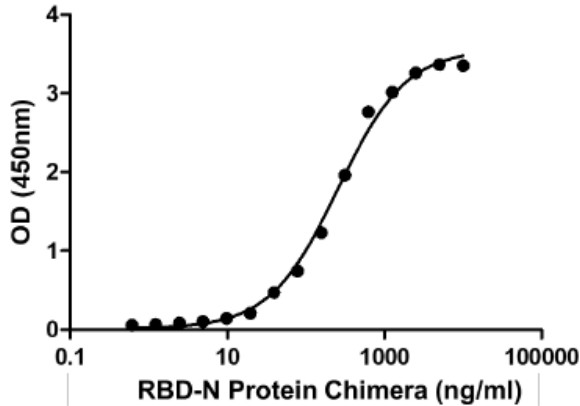
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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.comPlease 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

**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: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com