



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

Anti-Nucleocapsid Antibody (SARS-CoV-2)

Human, Recombinant, Monoclonal

Catalog #: 100861

Lot #: 200820

Conc: 0.20 mg/ml

Host Species | Isotype:

Human | Human IgG1

Formulated in: PBS and 0.02% sodium azide

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

References:

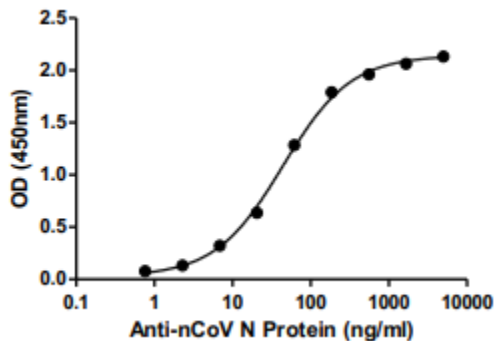
1. Surjit, M., *et al.*, *J. Biol. Chem.* 2006; **281**: 10669-10681.
2. Zeng, W., *et al.*, *Biochem. Biophys. Res. Commun.* 2020; **527(3)**: 618-623.

Description: Recombinant human monoclonal antibody recognizing the SARS-CoV-2 Nucleocapsid (N) protein.

Purification: Purified using protein A resin.

Applications: Useful in ELISA and Western blotting assays (1:1000) for detecting Nucleocapsid (N) protein.

Quality Assurance



Binding of Anti-Nucleocapsid Antibody (SARS-CoV-2) to Nucleocapsid Protein, Avi-His-tag (SARS-CoV-2) (BPS Bioscience #100778)

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**

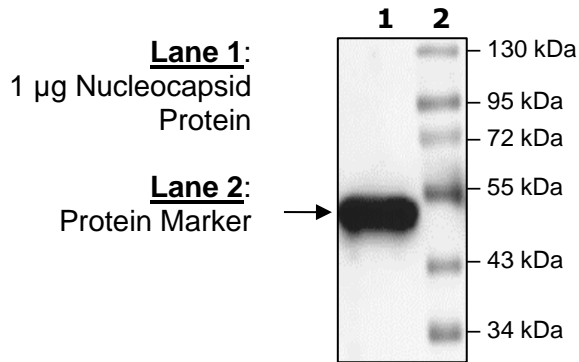
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Western Blot (1:1000)



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