



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

## Data Sheet

### Anti-SARS-CoV-2 Nucleocapsid Antibody

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.

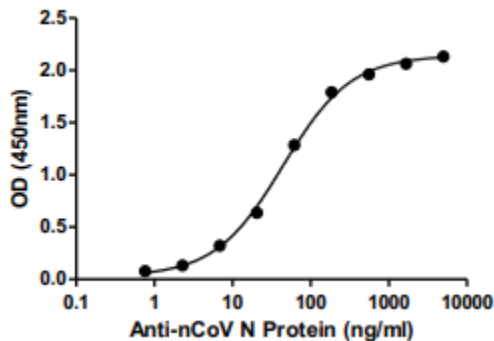
**Description:** Recombinant human monoclonal antibody recognizing the SARS-CoV-2 Nucleocapsid (N) protein.

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

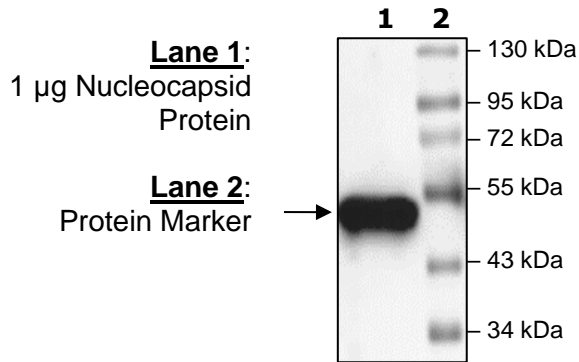
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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