

## Data Sheet

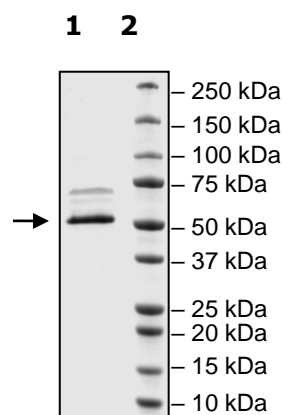
### **Nucleocapsid Protein, Avi-His-tag, Biotin-Labeled (SARS-CoV-2)**

SARS-CoV-2, Recombinant, C-terminal  
Avi-His-tag, Biotin-labeled**Catalog #:** 100779**Lot #:** 201008**Conc.:** 1.65 mg/ml**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3  
mM DTT.**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . *Avoid  
freeze/thaw cycles.***References:**

1. Wu, A. *et al.* Cell Host & Microbe.  
**27(3):** 325-328 (2020).
2. Benvenuto, D., *et al.* Medical  
Virology. **92(4):** 455-459 (2020).

**Description:** SARS-CoV-2 nucleocapsid protein, also known as COVID-19 nucleocapsid and SARS-CoV-2 N protein, Genbank Accession No.: YP\_009724397.2, a.a. 1-419(end), with C-terminal Avi-His-tag, expressed in a HEK293 cell expression system. MW= 48 kDa. This protein runs at a higher M.W. by SDS-PAGE due to glycosylation. This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be  $\geq 90\%$ 

## **Quality Assurance**

**4-20% SDS-PAGE  
Coomassie staining****Lane 1:  
Nucleocapsid****Lane 2:  
Protein Marker****Purity:**  $\geq 90\%$ **MW:** 48 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](mailto:info@bpsbioscience.com)Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)