

## Data Sheet

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

SARS-CoV-2, Recombinant, C-terminal  
Avi-His-tag

**Catalog #:** 100778

**Lot #:** 200601-1

**Conc.:** 1.07 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, and 20% glycerol.

**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.

## **Quality Assurance**

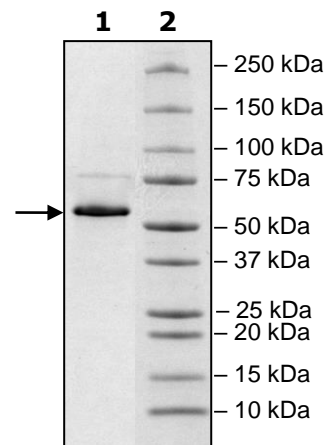
### 4-20% SDS-PAGE Coomassie staining

Lane 1:  
2 $\mu\text{g}$  Nucleocapsid

Lane 2:  
Protein Marker

MW: 48 kDa +glycans

Purity: >90%



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)