

## SARS-CoV-2 Nucleocapsid Protein

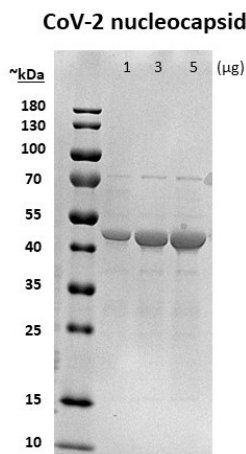
Cat. # CV2002

**Background:** Within the last two decades, SARS and MERS coronaviruses emerged as global health concerns causing severe acute respiratory syndromes. In December 2019, a novel coronavirus (SARS-CoV-2) was identified in Wuhan, Hubei province in China (1-3). The SARS-CoV-2 genome encodes several structural proteins including the nucleocapsid protein (Nucleoprotein; Protein N), which plays a role in virion assembly through its interaction with the viral genome and the membrane protein. This key protein packages the positive strand viral RNA into a helical ribonucleocapsid. The nucleocapsid represents a valuable tool for diagnostic and vaccine production purposes.

**Alternate names:** Nucleoprotein; Protein N, NC

### Product Information

<b>Molecular Weight:</b>	46 kDa (residues 1-419)
<b>Quantity:</b>	100 µg
<b>Physical State:</b>	Liquid
<b>Species:</b>	SARS-CoV-2
<b>Tag:</b>	None
<b>Activity:</b>	~99%
<b>Purity:</b>	~99%
<b>Storage:</b>	-80° C. Avoid repeated freeze/thaw cycles.



### References

1. Wang D, Hu B, Hu C, et al. Clinical Characteristics of 138 Hospitalized Patients With 2019 Novel Coronavirus–Infected Pneumonia in Wuhan, China. *JAMA*. 2020;323(11):1061.
2. Zhou P, Yang X-L, Wang X-G, et al. A pneumonia outbreak associated with a new coronavirus of probable bat origin. *Nature*. 2020;579(7798):270–273.
3. Zhu N, Zhang D, Wang W, et al. A Novel Coronavirus from Patients with Pneumonia in China, 2019. *N. Engl. J. Med.* 2020;382(8):727–733.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2009 LifeSensors, Inc. All Rights Reserved