

# SARS Nucleocapsid Protein Antibody (Protein N, Nucleoprotein)

Cat. # CV4001

---

## 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. The SARS-CoV 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 is a highly conserved and immunogenic viral protein thus representing a valuable tool for diagnostic and vaccine production purposes. The SARS Nucleocapsid Protein Antibody recognizes and binds to both SARS-CoV and SARS-CoV-2 nucleocapsid proteins.

**Target Molecular Weight:** 46 kDa

---

## Product Information

**Description:** Rabbit, polyclonal antibody to SARS Nucleocapsid Protein

**Species Cross Reactivity:** SARS-CoV, SARS-CoV-2

**Source:** Rabbit

**Applications:** WB, ELISA

**Recommended Antibody Dilutions:**

Western Blotting: Robust detection of 10 ng of recombinant protein was possible when antibody was used at a final concentration of 1 µg/mL

---

## Storage/Purification

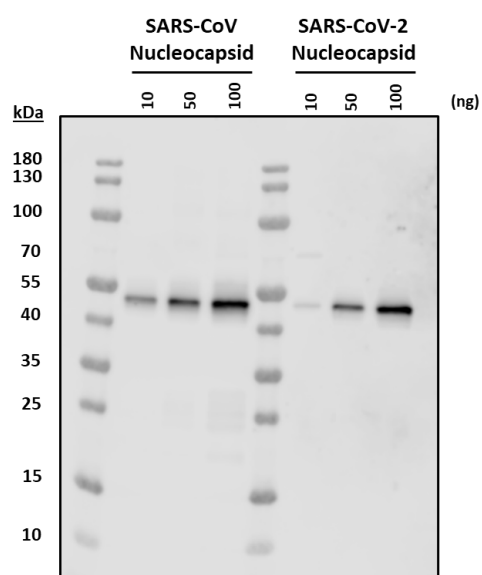
Polyclonal antibodies are produced by repeatedly immunizing rabbits with purified recombinant full-length protein. Antibodies are purified from monospecific antiserum by protein A affinity purification.

**Storage:** Supplied in 0.02 M Potassium Phosphate, 0.15 M Sodium Chloride pH 7.2.

Store at -20°C.

*Avoid Freeze/Thaw Cycles.*

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2007 LifeSensors, Inc. All Rights Reserved



## 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 © 2007 LifeSensors, Inc. All Rights Reserved