

---

## SARS-CoV-2 N Protein Monoclonal Antibody

(Catalog # A73663)

---

### Background

Coronaviruses are enveloped viruses with a positive-sense RNA genome and with a nucleocapsid of helical symmetry. Coronavirus nucleoproteins localize to the cytoplasm and the nucleolus, a subnuclear structure, in both virus-infected primary cells and in cells transfected with plasmids that express N protein. Coronavirus N protein is required for coronavirus RNA synthesis, and has RNA chaperone activity that may be involved in template switch. Nucleocapsid protein is a most abundant protein of coronavirus. During virion assembly, N protein binds to viral RNA and leads to formation of the helical nucleocapsid. Nucleocapsid protein is a highly immunogenic phosphoprotein also implicated in viral genome replication and in modulating cell signaling pathways. Because of the conservation of N protein sequence and its strong immunogenicity, the N protein of coronavirus is chosen as a diagnostic tool.

### Description

SARS-CoV-2 N Protein Monoclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Buffer: PBS with 0.02% sodium azide, pH7.3.

### Specificity

2019-nCoV

### Isotype

IgG

### Uniprot ID

P59595

### Purification

Affinity Purification

### Immunogen

Recombinant fusion protein of SARS-COV-2 Nucleoprotein.

### Storage

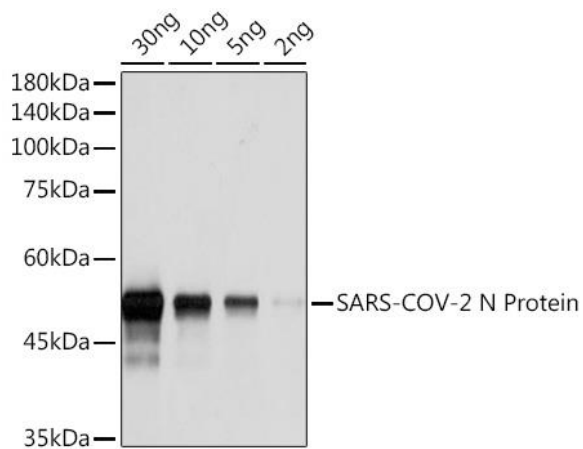
Shipped at 4°C. Store at 4°C. Avoid freeze / thaw cycles.

### Alternative Names

Anti-coronavirus NP Antibody, Anti-coronavirus Nucleocapsid Antibody, Anti-coronavirus Nucleoprotein Antibody, Anti-cov np Antibody, Anti-ncov NP Antibody, Anti-novel coronavirus NP Antibody, Anti-novel coronavirus Nucleocapsid Antibody, Anti-novel coronavirus Nucleoprotein Antibody, Anti-NP Antibody, Anti-Nucleocapsid Antibody, Anti-Nucleoprotein Antibody

### Application

WB, ELISA, FCM, IHC, IF, IP; Recommended dilution: WB, 1:1000 - 1:5000; ELISA, 1:1000 - 1:5000; FCM, 1:50 - 1:200; IHC, 1:50 - 1:200; IF, 1:50 - 1:200; IP, 1:50 - 1:200



Western blot analysis of SARS-CoV-2 N Protein Monoclonal Antibody at 1:1000 dilution.

Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 5s.