

---

## SARS-CoV-2 Neutralizing Recombinant Monoclonal Antibody [DA0002]

(Catalog # A73683)

---

### Description

SARS-CoV-2 Neutralizing Recombinant Monoclonal Antibody is isolated from a SARS-CoV-2 infected patient and is recombinantly produced from human 293 cells (HEK293). This antibody recognizes the SARS-CoV-2 Spike Protein RBD domain and inhibits the interaction between SARS-CoV-2 RBD and ACE2. Clone: DA0002

### Formulation

Liquid. Lyophilized from a 0.22 µm filtered solution of PBS, pH 7.2, with 10% Trehalose.

### Host

SARS-CoV-2

### Source

Human

### Specificity

SARS-CoV-2

### Isotype

IgG1 (Human)

### Purification

Affinity Purification; >95% by SDS-PAGE

### Endotoxin

< 0.1 EU/µg of the protein by LAL method.

### Storage

Store the lyophilized protein at -20°C to -80 °C for long term. Avoid repeated freeze/thaw cycles.

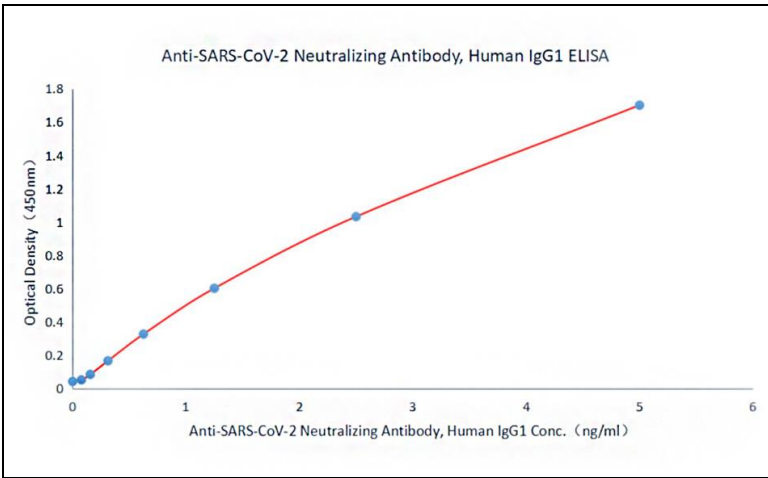
### Alternative Names

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; SARS-CoV-2; 2019-nCoV

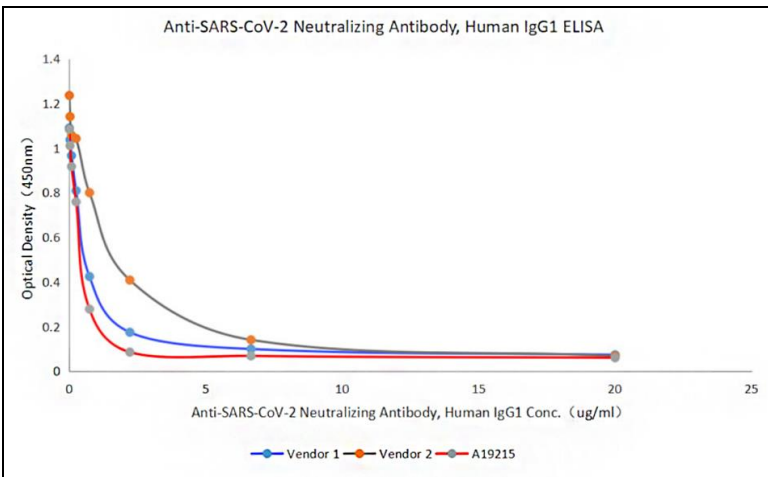
### Application

ELISA: Recommended dilution: ELISA 1:250 – 1:1000.

---



Immobilized SARS-CoV-2 S1 protein, His Tag (Cat# E80004-1) at 1 µg/mL (100µL/well) can bind Anti-SARS-CoV-2 Neutralizing Antibody, Human IgG1 with a linear range of 0.078-5 ng/mL.



Serial dilutions of Anti-SARS-CoV-2 Neutralizing Antibody, Human IgG1 was detected with a half maximal inhibitory concentration (IC50) of 0.496ug/mL.