

SARS-CoV-2 Nucleocapsid IgG ELISA Kit (RUO)

SARS-CoV-2 Nucleocapsid IgG ELISA kit enabling detection of human SARS-CoV-2 IgG in serum and plasma.

The SARS-CoV-2 Nucleocapsid IgG ELISA Kit is designed for measuring the amount of SARS-CoV-2 nucleocapsid IgG in serum and plasma samples with only 30 minutes in incubation time. Absorbance is read at 450 nm. Measure up to 86 samples with 1 kit.

Application Note: [Qualitative Measurement of SARS-CoV-2 Spike and Nucleocapsid IgG Antibodies in Serum Samples](#)

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

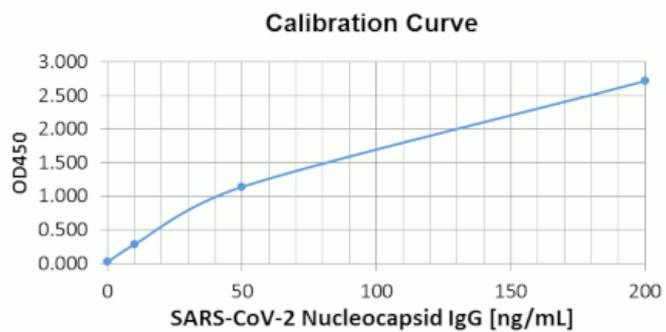
ENZ-KIT193-0001

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Sensitive detection of IgG antibodies against SARS-CoV-2 nucleocapsid protein
- Rapid qualitative or semi-quantitative results in 30 minutes
- High-throughput immunoassay measures up to 86 samples in singlicate or 43 in duplicate.



Typical Calibration Curve for SARS-CoV-2
Nucleocapsid IgG ELISA Kit

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Coronavirus (COVID-19) IgG
Application	Colorimetric detection, ELISA
Application Notes	For the semi-quantitative determination of human SARS-CoV-2 IgG in serum.
Assay Time	30 minutes
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Contents	SARS-CoV-2 Nucleocapsid Coated Microplate SARS-CoV-2 Nucleocapsid IgG Calibrator Pack SARS-CoV-2 Nucleocapsid Conjugate SARS-CoV-2 Nucleocapsid Sample Diluent Wash Buffer Concentrate (20X) TMB Substrate Stop Solution 2 SARS-CoV-2 Nucleocapsid Assay Layout Sheet
Species Reactivity	Human
Wavelength	450 nm

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com