

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

Highly sensitive SARS-CoV-2 Spike IgG ELISA kit enabling detection of human SARS-CoV-2 IgG in serum.

The SARS-CoV-2 Spike IgG ELISA kit is a colorimetric immunoassay kit with results in under 2 hours. Absorbance is read at 450 nm. Measure up to 86 samples with 1 kit.

Citations: 1

[View Online »](#)

Ordering Information

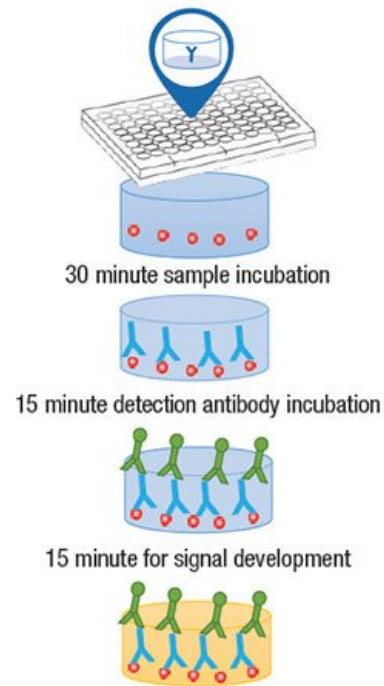
[Order Online »](#)

ENZ-KIT190-0001	96 wells
-----------------	----------

Manuals, SDS & CofA

[View Online »](#)

- High-sensitivity ELISA enables detection of low amounts of SARS-CoV-2 Spike S1-RBD IgG
- Rapid qualitative results in under 2 hours
- High-throughput immunoassay measures up to 86 samples per kit



Schematic diagram of the assay protocol for the SARS-CoV-2 IgG ELISA Kit.

Handling & Storage

Long Term Storage +4°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 qualitative determination of human SARS-CoV-2 IgG in serum.

Assay Time <2 hours

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents SARS-CoV-2 Antigen Coated Microplate
High Positive Control
Low Positive Control
Negative Control
Sample Diluent
Wash Buffer Concentrate
Stop Solution
TMB Substrate
HRP Conjugate
Plate Sealer

Species Reactivity Human

Wavelength 450 nm

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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