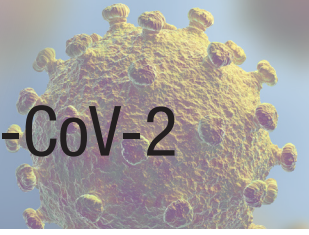




# Antibody Detection for SARS-CoV-2



## SARS-CoV-2 IgG ELISA Kit

Our Enzo SARS-CoV-2 IgG ELISA Kit is an Enzyme-Linked Immunosorbent Assay (ELISA) designed for the qualitative detection of IgG antibodies specific to SARS-CoV-2 in human serum samples issued for use under Emergency Use Authorization (EUA) as an aid in identifying individuals with an adaptive immune response to SARS-CoV-2.

Formatted in a 96-well microplate (12 x 8 well strips), the assay is easily adapted for either automated open platforms or manual workflows.

Designed for accurate and sensitive detection, the indirect ELISA has a two-step binding process involving a specific SARS-CoV-2 antigen and a HRP conjugated anti-human IgG secondary antibody.

### HIGH PERFORMANCE

Clinical Specificity\*: 96.5%

Clinical Sensitivity\*: 100%

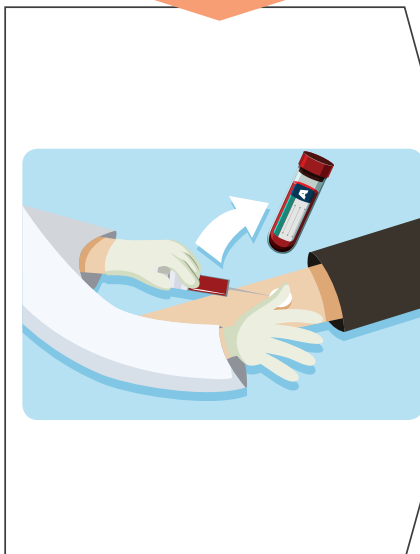
Time to results: < 2 hrs

\*In a clinical study with a RT-PCR comparator

## SARS-CoV-2 IgG ELISA Kit (ENZ-KIT170)

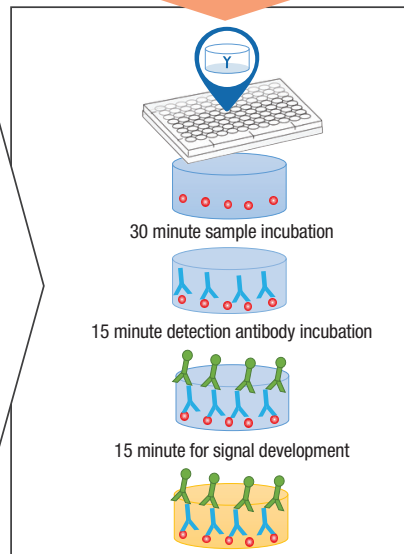
TIME TO RESULT: < 2 HOURS

### Sample Collection



Serum Sample

### Antibody Detection



Indirect ELISA

### Qualitative Analysis



Positive or Negative IgG

## SARS-CoV-2 IgG ELISA Kit (ENZ-KIT170)

COMPONENTS	QUANTITY	STORAGE
SARS-CoV-2 Antigen Coated Microplate	1 x 96-well plate	4°C
High Positive Control	1 x 500 µL	4°C
Low Positive Control	1 x 500 µL	4°C
Sample Diluent	1 x 10 mL	4°C
Wash Buffer Concentrate (20X)	1 x 50 mL	4°C
Stop Solution	1 x 10 mL	4°C
TMB Substrate	1 x 10 mL	4°C
HRP Conjugate	1 x 10 mL	4°C
Plate Sealer	2 each	–

### RELATED PRODUCTS

Product Name	Product Number
AMPIXTRACT™ SARS-CoV-2 Extraction Kit	ENZ-GEN216
AMPIPROBE® SARS-CoV-2 RT-PCR Assay Kit	ENZ-GEN215
AMPIPROBE® SARS-CoV-2 Controls	ENZ-GEN218
IFN-γ (human) ELISA Kit	ADI-900-136
IL-1β (human) ELISA Kit	ADI-900-130A
IL-6 (human) High Sensitivity ELISA Kit	ENZ-KIT178
IL-6 (human) Sensitivity Controls	ENZ-KIT178-CTL
TNF-α (human) ELISA Kit, 1 Pack	ADI-900-099
TNF-α (human) ELISA Kit, 5 Pack	ADI-901-099

**Global Headquarters**  
**ENZO LIFE SCIENCES, INC.**  
 Ph: 800.942.0430  
 info-usa@enzolifesciences.com

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

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

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

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

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

For local distributors and detailed product information visit us online: [www.enzolifesciences.com/coronavirus](http://www.enzolifesciences.com/coronavirus)

For *In Vitro* Diagnostic Use. For Emergency Use Authorization Only.