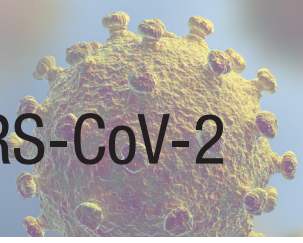




Antibody Detection for SARS-CoV-2



SARS-CoV-2 Spike Protein IgG ELISA Kit

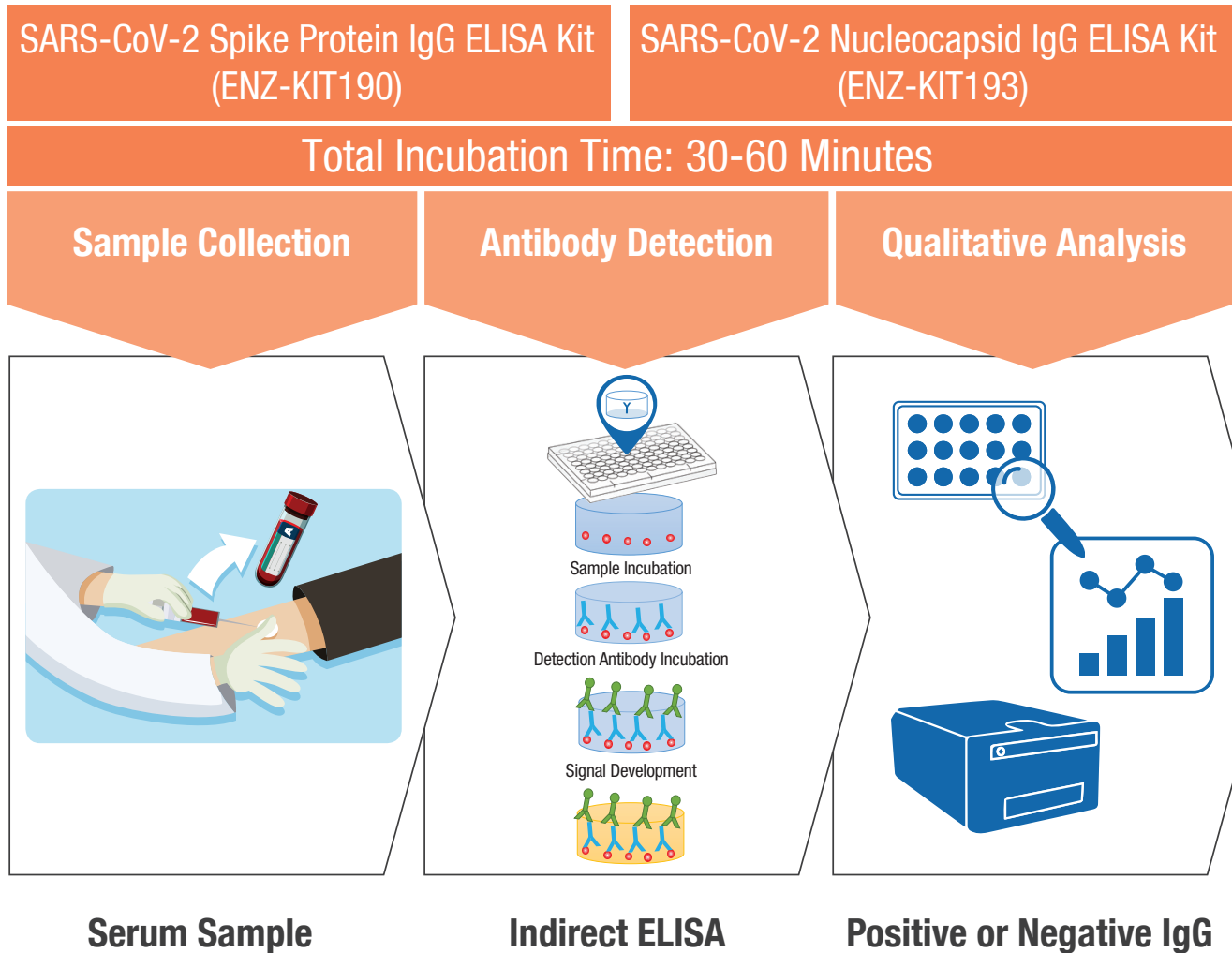
RUO

SARS-CoV-2 Nucleocapsid IgG ELISA Kit

RUO

Our Enzo SARS-CoV-2 IgG ELISA Kit featuring a SARS-CoV-2 Spike Protein Coated Microplate and SARS-CoV-2 Nucleocapsid IgG ELISA Kit featuring a SARS-CoV-2 Nucleocapsid Protein Coated Microplate are Enzyme-Linked Immunosorbent Assays (ELISAs) designed for the qualitative detection of IgG antibodies specific to SARS-CoV-2 in human serum samples as an aid in identifying individuals with an adaptive immune response to SARS-CoV-2 indicating recent or prior infection. Formatted in a 96-well microplate, 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.



Serum Sample

Indirect ELISA

Positive or Negative IgG

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

ENZ-KIT190

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

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

ENZ-KIT193

COMPONENTS	QUANTITY	STORAGE
SARS-CoV-2 Nucleocapsid Coated Microplate	1 x 96-well plate	Long Term: -20°C Short Term (1 month): 4°C
Calibrator Pack	Calibrators 1-4	
Conjugate	1 x 10 mL	
Sample Diluent	1 x 100 mL	
Wash Buffer Concentrate (20X)	1 x 30 mL	
Stop Solution	1 x 10 mL	
TMB Substrate	1 x 10 mL	

RELATED PRODUCTS

Product Name	Product Number
AMPIXTRACT™ SARS-CoV-2 Extraction Kit (RUO)	ENZ-GEN225
AMPIPROBE® SARS-CoV-2 Assay Kit (RUO)	ENZ-GEN230
AMPIPROBE® SARS-CoV-2 Controls (RUO)	ENZ-GEN231

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 detailed product information visit us online: www.enzolifesciences.com/coronavirus

For Research Use Only.

All product names, logos, and brands are the property of their respective owners.