

# Coronavirus (COVID-19, Spike) polyclonal antibody

This antibody is covered by our [Worry-Free Guarantee](#).

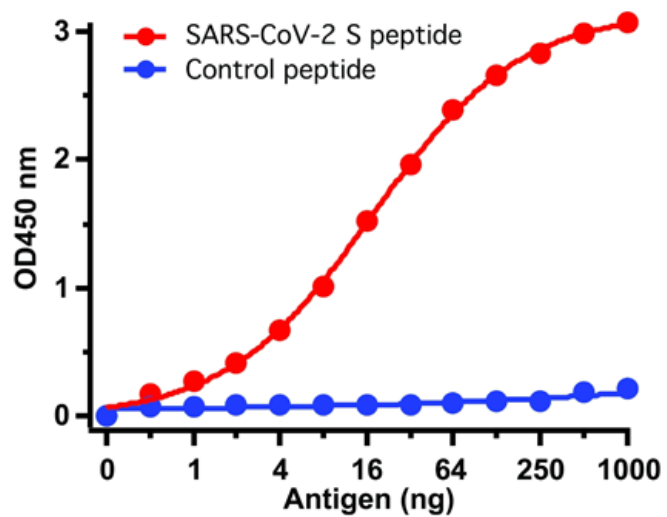
## Ordering Information

[Order Online »](#)

ENZ-ABS827-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



ELISA using SARS-CoV-2 (COVID-19, 2019-nCoV)  
Spike Antibody

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** SARS-CoV-2 Spike

**Application** ELISA

**Formulation** Liquid. In PBS containing 0.02% sodium azide.

**Host** Rabbit

**Immunogen** Peptide corresponding to 20 amino acids near the carboxy terminus of SARS-CoV-2 (COVID-19, 2019-nCoV) Spike glycoprotein. The immunogen is located within the last 50 amino acids of SARS-CoV-2 (COVID-19, 2019-nCoV) spike protein.

**Isotype** IgG

**Purity Detail** Affinity purified.

**Species Reactivity** Coronavirus

**Worry-free Guarantee** This antibody is covered by our [Worry-Free Guarantee](#).

**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-  
be@enzolifesciences.com](mailto:info-be@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)