## Coronavirus (SARS-CoV Spike) polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

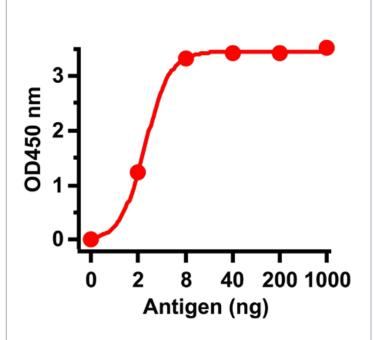
Order Online »

ENZ-ABS825-0100

100µg

Manuals, SDS & CofA

**View Online »** 



A sandwich ELISA was performed using immunogen as coating antigen and the SARS-CoV Spike antibody as the capture antibody.

## **Handling & Storage**

**Handling** Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

**Application** ELISA

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

**Host** Rabbit

**Immunogen** Peptide corresponding to 17 amino acids near the center

of SARS-CoV Spike glycoprotein. The immunogen is located within amino acids 550-600 of SARS-CoV Spike.

**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

