

Coronavirus (SARS-CoV S1) antigen (aa 1-53, 90-115, 171-205), (recombinant) (His-tag)

Ordering Information

[Order Online »](#)

ENZ-PRT299-0050	50µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

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 Coronavirus spike protein

Formulation Liquid. In 10 mM PBS, pH 7.2.

Purity ≥90%

Purity Detail Purified.

Source Produced in *E. coli*. Coronavirus (SARS-CoV) spike protein (aa 1-53, 90-115, 171-205) is fused to a His-tag.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

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

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

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