

Coronavirus (SARS-CoV S1) antigen (aa 12-53, 90-115, 171-203), (recombinant)

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 25 mM Tris-HCl with detergents and 50% glycerol.
MW	38 kDa
Purity	≥95%
Purity Detail	Purified.
Source	Produced in <i>E. coli</i> . Coronavirus (SARS-CoV) spike protein (aa 12-53, 90-115, 171-203).

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