

Coronavirus (SARS-CoV S2) antigen (aa 408-470, 540-573), (recombinant) (His-tag)

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 <i>E. coli</i> . Coronavirus (SARS-CoV) spike protein (aa 408-470, 540-573) 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