

# Coronavirus (MERS-CoV S1) antigen, (recombinant) (His-tag)

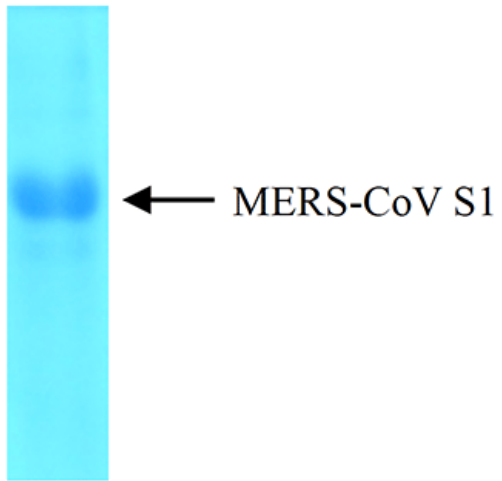
## Ordering Information

[Order Online »](#)

ENZ-PRT309-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Coomassie-stained SDS-PAGE showing purified recombinant MERS-CoV Spike (S1) protein

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -80°C

**Shipping** Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Formulation** Liquid. In PBS, 25mM Arginine and 0.05% Sodium Azide

**Purity** ≥ 95%

**Purity Detail** Purified.

**Source** Produced in *E. coli*. Recombinant protein (aa 56-295).

**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)