

# Zika Virus -NS1 (Duo pack)

Recombinant Zika Virus NS1 protein, with C-terminal 6x His tags, produced in mammalian HEK293 cells. Includes one vial each of Zika NS1 from Suriname and Uganda strains

**Manuals, SDS & CofA**

[View Online »](#)

## Handling & Storage

Long Term Storage -80°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	ZIKV-NS1
Formulation	Liquid. In DPBS pH 7.4.
Purity	≥85% (SDS-PAGE)
Purity Detail	Purified protein.
Source	Produced in mammalian HEK293 cells. Recombinant Zika Virus NS1 protein, with C-terminal 6x His tags.

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)