

Zika Virus -NS1 (Suriname strain)

Recombinant Zika Virus (Suriname strain) NS1 protein, incorporating a C-terminal 6x His tag produced in mammalian HEK293 cells.

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	≥90% (SDS-PAGE)
Purity Detail	Purified protein.
Source	Produced in mammalian HEK293 cells. Recombinant Zika Virus (Suriname strain) NS1 protein, incorporating a C-terminal 6x 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