

Zika Virus Envelope protein (Suriname strain)

Zika virus envelope protein recombinant antigen containing a C-terminal His-tag

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Short Term Storage +4°C

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Formulation Liquid. In 20mM NaH₂PO₄ pH 7.8, 210mM NaCl.

Purity ≥95% (SDS-PAGE)

Purity Detail Purified protein.

Source Produced in HEK293 mammalian cells. Recombinant Zika virus envelope protein with C-terminal 6-His tag.

UniProt ID Q91KX7

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