

Zika Virus VLP

Recombinant Zika virus-like particles consisting of Envelope, pre-Membrane, and Membrane protein produced in mammalian HEK293 cells.

Ordering Information

[Order Online »](#)

ENZ-PRT278-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	ZIKV-VLP
Appearance	Light-brown opaque liquid.
Formulation	Liquid. In DPBS pH 7.4, 20% Sucrose. May contain traces of tartrate.
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified protein.
Source	Produced in mammalian HEK293 cells. Recombinant Zika virus-like particles consisting of envelope, pre-membrane, and membrane protein.

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