

Dengue NS1 ST3, (recombinant)

Dengue virus NS1 protein serotype 3 presented in its native folded state with all post-translational modifications to deliver optimal antigenicity.

Ordering Information

[Order Online »](#)

ENZ-PRT106-0100	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 DENV1-NS1

Appearance Clear liquid.

Formulation Liquid. In DPBS pH 7.4.

Purity ≥95% (SDS-PAGE)

Purity Detail Purified protein.

Source Produced in HEK 293 cells. Dengue Viral NS1 Serotype 3 (DENV1-NS1) from strain Sri Lanka D3/H/IMTSSA-SRI/2000/1266.

Technical Info / Product Notes We offer all four Dengue Virus serotypes.

UniProt ID Q6YMS4



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