

Dengue NS1 ST1, (recombinant)

Dengue virus NS1 protein serotype 1 presented in its native folded state with all post-translational modifications to deliver optimal antigenicity.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-PRT104-0100	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 1 (DENV1-NS1) from strain Nauru/Western Pacific/1974.
Technical Info / Product Notes	We offer all four Dengue Virus serotypes.
UniProt ID	P17763

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