

# Dengue NS1 ST2, (recombinant)

Dengue virus NS1 protein serotype 2 presented in its native folded state with all post-translational modifications to deliver optimal antigenicity.

Citations: 2

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-PRT105-0100	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	DENV2-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 2 (DENV2-NS1) from strain Thailand/16681/84.
Technical Info / Product Notes	We offer all four Dengue Virus serotypes.
UniProt ID	P29990

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)