

# Zika Virus NS1 monoclonal antibody (D11)

This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Application	ELISA, IF
Clone	D11
Crossreactivity	It demonstrates negligible cross-reactivity with NS1 proteins from Dengue virus (all serotypes), Japanese Encephalitis Virus and Yellow Fever Virus. A small amount of cross-reactivity has been observed with the NS1 protein from West Nile Virus in a direct ELISA.
Formulation	Liquid. In PBS, pH 7.4.
Host	Mouse
Isotype	IgG1
Purity Detail	Protein G affinity purified.
Source	Purified from hybridoma tissue culture supernatant.
Specificity	Detects NS1 from both the Uganda and Suriname strains.
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

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-fr@enzolifesciences.com](mailto:info-fr@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)