

# Tick-borne Encephalitis Virus NS1 monoclonal antibody (M834)

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

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Application	ELISA, WB
Clone	M834
Crossreactivity	In ELISA assays, the antibody does not cross-react with other flavivirus NS1 proteins tested including Yellow Fever Virus (YFV), Japanese Encephalitis Virus (JEV), Powassan Virus (PV), Zika virus or DENV (a mix of DENV1, DENV2, DENV3, DENV4). In Western Blotting, the antibody predominantly recognises the dimeric form of TBEV NS1. The antibody also cross reacts with PV NS1, JEV NS1 and DENV NS1 but not with YFV or ZIKV NS1.
Formulation	Liquid. In PBS, pH 7.2, containing 0.05% sodium azide.
Host	Mouse
Isotype	IgG1
Source	Purified from hybridoma tissue culture supernatant.
Specificity	Recognizes Tick-borne Encephalitis Virus (TBEV) NS1 protein.
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-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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

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