Tick-borne Encephalitis Virus NS1 monoclonal antibody (M834)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS837-0100

100µg

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

lgG1

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 Worry-Free Guarantee

.

