Mu opioid conformation specific polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS695-0100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA, Flow Cytometry, IHC, IP, WB

Application NotesSuitable for cell-based and membrane-based ELISA.

Predicted MW of ~44kDa.

Formulation Liquid. In PBS containing 50% glycerol and 0.01% sodium

azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to the N-terminus of the

Mu opioid receptor (SDPLAPASWSPAPGSWL).

Purity Detail Affinity purified.

Recommendation Dilutions/Conditions Conformation ELISA (1:500)TITER by direct ELISA

(1:1000)Immunohistology (1:1000)Western Blot (1:750)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be

determined individually for each application.

Species Reactivity Human, Mouse, Rat

UniProt ID P33535 / P35372

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com