Netrin-1 monoclonal antibody (Nora-1)

This antibody is covered by our Worry-Free Guarantee.

Citations: 6

View Online »

Ordering Information

Order Online »

ALX-804-838-C100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability Stable for at least 1 year after receipt when stored at -20°C.

Handling Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA, IHC, IP, WB

Clone Nora-1

Formulation Liquid. In PBS containing 0.02% sodium azide.

Host Mouse

Immunogen Recombinant human, mouse and chicken netrin-1 (aa 28-

264).

lgG2a

Positive Control Prod. No. ALX-522-100

Recommendation Dilutions/Conditions Immunoprecipitation (1:100)Western Blot

(1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be

determined individually for each application.

Source Purified from concentrated hybridoma tissue culture

supernatant.

Species Reactivity Chicken, Human, Mouse

UniProt ID 095631

Last modified: May 29, 2024



info-

eu@enzolifesciences.com