TWEAK (human) monoclonal antibody (Mira-2)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ALX-804-842-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 reconstitution, 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

Alternative Name Apo-3L, TNFSF 12

Application ELISA, Flow Cytometry, IP

Application Notes Immunoprecipitation (1:100)

Optimal conditions must be determined individually for

each application.

Clone Mira-2

Formulation Liquid. In PBS containing 0.02% sodium azide.

Host Mouse

Immunogen Recombinant human TWEAK extracellular domain (aa

106-249) (Prod. No. ALX-522-021).

lgG1

Prod. No. ALX-522-021

Purity Detail Protein G-affinity purified.

Recommendation Dilutions/Conditions Immunoprecipitation (1:100)Suggested dilutions/conditions

may not be available for all applications. Optimal conditions

must be determined individually for each application.

Purified from concentrated hybridoma tissue culture Source

supernatant.

Species Reactivity Human

UniProt ID O43508

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

Last modified: May 29, 2024



info-

uk@enzolifesciences.com