TRAIL (human) monoclonal antibody (III6F)

This antibody is covered by our Worry-Free Guarantee.

Citations: 5

View Online »

Ordering Information

Order Online »

ALX-804-326-C100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability Stable for at least 1 year after receipt when stored as recommended.

Handling Avoid freeze/thaw cycles.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Apo-2L, TNFSF 10, CD253

Application ELISA, Flow Cytometry, IHC (FS), IP, WB

Clone III6F

Formulation Liquid. In PBS containing 0.02% sodium azide.

GenBank ID U57059

Gene/Protein Identifier 8743 (Entrez GenelD)

Host Mouse

Immunogen Recombinant human soluble TRAIL.

Isotype IgG2b

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human

UniProt ID P50591

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

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com