DCC (human) monoclonal antibody (Axel-1)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ALX-804-853-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.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Deleted in colorectal cancer protein

Application ELISA, Flow Cytometry, FUNC

Clone Axel-1

Formulation Liquid. In PBS containing 0.02% sodium azide.

Host Mouse

Immunogen Recombinant human DCC fibronectin domain 5 (Prod. No.

ALX-522-109).

lgG2a λ

Purity Detail Protein G-affinity purified.

Source Purified from concentrated hybridoma tissue culture

supernatant.

Species Reactivity Human

UniProt ID P43146

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

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com