CD3 (human) monoclonal antibody (MEM-57)

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

Ordering Information

Order Online »

ALX-805-057-C100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability Stable for 6 months when stored at +4°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 Flow Cytometry, FUNC, IP

Application Notes Functional Application: activates T cells *in vitro*; highly

effective immunosuppressive agent.

Clone MEM-57

Formulation Liquid. In PBS, pH 7.4, containing 15mM sodium azide.

Host Mouse

Immunogen Human thymocytes and T lymphocytes.

lgG2a

Purity Detail Protein A-affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (2-5µg/ml))Suggested dilutions/conditions

may not be available for all applications. Optimal conditions

must be determined individually for each application.

Species Reactivity Human

Specificity Recognizes the ε and δ chain of human CD3.

UniProt ID P09693 (γ chain), P04234 (δ chain), P07766 (ε chain),

P20963 (ζ chain)

Last modified: May 29, 2024



info-

eu@enzolifesciences.com