

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 <i>in vitro</i> ; 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.
Isotype	IgG2a
Purity Detail	Protein A-affinity purified.
Recommendation	Flow Cytometry (2-5µg/ml))Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Dilutions/Conditions	
Species Reactivity	Human
Specificity	Recognizes the ε and δ chain of human CD3.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
P09696 (γ chain), P04234 (δ chain), P07766 (ε chain), P20963 (ζ chain)
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
P09696 (γ chain), P04234 (δ chain), P07766 (ε chain), P20963 (ζ chain)
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
P09696 (γ chain), P04234 (δ chain), P07766 (ε chain), P20963 (ζ chain)
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
P09696 (γ chain), P04234 (δ chain), P07766 (ε chain), P20963 (ζ chain)
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

Worry-free Guarantee

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

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.