

CD3 ϵ Recombinant monoclonal antibody (500A2) (Mouse IgG1 κ)

This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|---------------------------|---|
| Use/Stability | Stable at 4°C for up to 3 months. For longer storage, aliquot and store 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 | T-cell surface glycoprotein CD3 ϵ chain, T-cell surface antigen T3/Leu-4 ϵ chain, CD3e |
| Application | Flow Cytometry |
| Application Notes | Functional Application: Activation |
| Clone | 500A2 |
| Formulation | Liquid. In PBS containing 0.02% Proclin 300. |
| Host | Mouse |
| Immunogen | T cell receptor complexes derived from mouse C6VL-BS thymoma cells (C57BL/Ka) . |
| Isotype | IgG1 κ |
| Purity Detail | Protein A affinity purified. |
| Species Reactivity | Mouse |
| Specificity | Recognizes the 25-kDa ϵ chain of the T-cell receptor-associated CD3 complex expressed on thymocytes, mature T lymphocytes, and NK-T cells on many mouse strains. |

Technical Info / Product Notes

This is a chimeric mouse antibody, based on the original hamster IgG format, created for improved compatibility with existing reagents, assays and techniques.

UniProt ID

P22646

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[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)