

CD45 (human) monoclonal antibody (MEM-28)

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

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-805-0410	100µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Flow Cytometry, ICC, IHC (PS), IP, WB

Clone MEM-28

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

Host Mouse

Immunogen Human thymocytes and T lymphocytes.

Isotype IgG1

Purity Detail Protein A-affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (1µg/ml)Immunocytochemistry (10µg/ml, paraformaldehyde fixation can be used)Immunohistochemistry (paraffin sections, no pre-treatment of tissue sections is essential)Western Blot (1µg/ml; Positive control: JURKAT human leukemia T-cell lysate, Kg-1a human leukemia cell lysate. Sample preparation: buffer with laurylmaltoside, 2x non-reducing SDS.Application note: non-reducing conditions. SDS-PAGE (6% separating gel))Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human

Specificity Recognizes all isoforms of CD45.

UniProt ID P08575

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)