

# CD45 (human) monoclonal antibody (MEM-28)

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

Citations: 5

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-805-041-C100	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

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)