

CD18 (human) monoclonal antibody (MEM-48)

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

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-805-062-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, 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

Alternative Name	β2 integrin
Application	Flow Cytometry, FUNC, IHC (FS), IP, WB
Application Notes	Functional Application: induces high-affinity LFA-1 conformation in leukocytes
Clone	MEM-48
Crossreactivity	Cross-reacts with all three leukocyte integrins.
Formulation	Liquid. In PBS, pH 7.4, containing 15mM sodium azide.
Host	Mouse
Immunogen	Leukocytes from a patient suffering from a LGL-type leukemia.
Isotype	IgG1
Purity Detail	Protein A-affinity purified.

Recommendation Dilutions/Conditions

Immunohistochemistry (paraffin sections: recommended dilution 10µg/ml, positive tissue: spleen, microglia; frozen sections)Western Blot (Recommended dilution: 2µg/ml in buffer containing laurylmaltoside, 2x non-reducing SDS; SDS-PAGE (6% separating gel, 4% stacking gel)positive control: Kg-1a human leukemia cell lysate)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity

Human

Specificity

Recognizes residues 534-546 in cysteine-rich repeat 3 of CD18 (β2 integrin).

UniProt ID

P05107

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)