

CD18 (human) monoclonal antibody (MEM-148)

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

Citations: 3

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Alternative Name	β2 integrin
Application	Flow Cytometry, FUNC, IP, WB
Application Notes	Functional Application: induces homotypic cell aggregation and high-affinity LFA-1 conformation in resting leukocytes.
Clone	MEM-148
Crossreactivity	Does not cross-react with pig CD18.
Formulation	Liquid. In PBS, pH 7.4, containing 15 mM sodium azide.
Host	Mouse
Immunogen	Human peripheral blood mononuclear cells.
Isotype	IgG1
Purity Detail	Protein A-affinity purified.
Recommendation Dilutions/Conditions	Western Blot (non-reducing)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
Specificity	Recognizes free CD18 (β2 integrin) not associated with CD11.

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)