

β 2-Microglobulin (human) monoclonal antibody (B2M-01)

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-4820	100 μ g
--------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Do not freeze.
Short Term Storage	+4°C
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	ELISA, Flow Cytometry, IP, WB
Clone	B2M-01
Crossreactivity	Does not cross-react with mouse, bovine, rabbit, dog or chicken β 2-microglobulin.
Formulation	Liquid. In PBS, pH 7.4, containing 15mM sodium azide.
Host	Mouse
Immunogen	Purified human β 2-microglobulin.
Isotype	IgG2a
Recommendation Dilutions/Conditions	Flow Cytometry (1 μ g/ml)Immunohistochemistry (paraffin sections, 10 μ g/ml)Western Blot (2-4 μ g/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from ascites by DEAE-chromatography and precipitation methods.
Species Reactivity	Human
Specificity	Recognizes human β 2-microglobulin associated with cell-surface MHC Class I molecules and other membrane antigens as well as soluble human β 2-microglobulin.

UniProt ID

P61769

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)