

β2-Microglobulin (human) monoclonal antibody (B2M-01)

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-482-C100	100μg
------------------	-------

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 Flow Cytometry (1 μ g/ml) Immunohistochemistry (paraffin sections, 10 μ g/ml) Western Blot
Dilutions/Conditions (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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +32 3 466 0420
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee