

β2-Microglobulin monoclonal antibody (34)

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

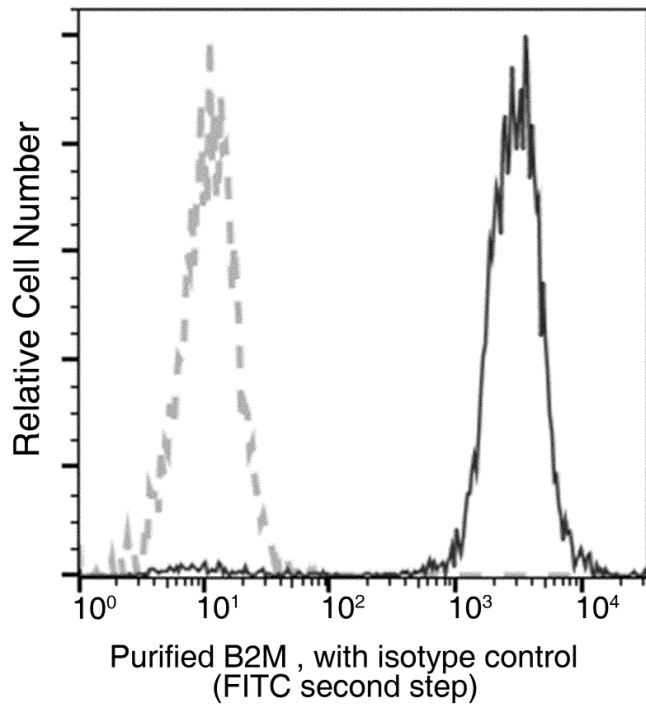
Ordering Information

[Order Online »](#)

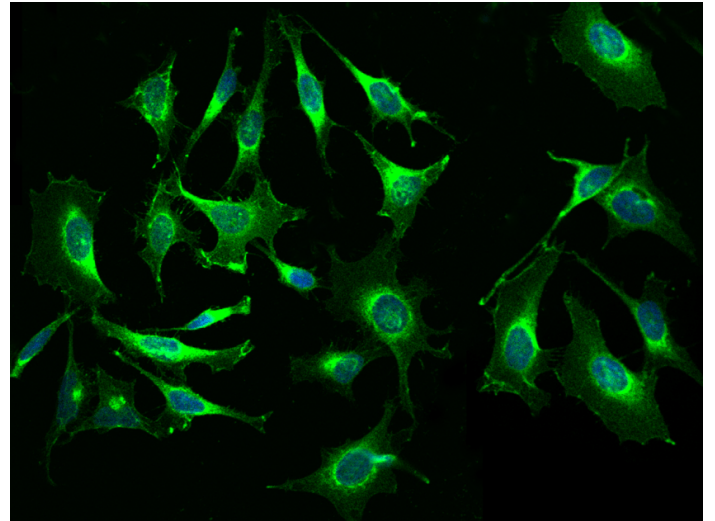
ENZ-ABS987-0100	100μl
-----------------	-------

Manuals, SDS & CofA

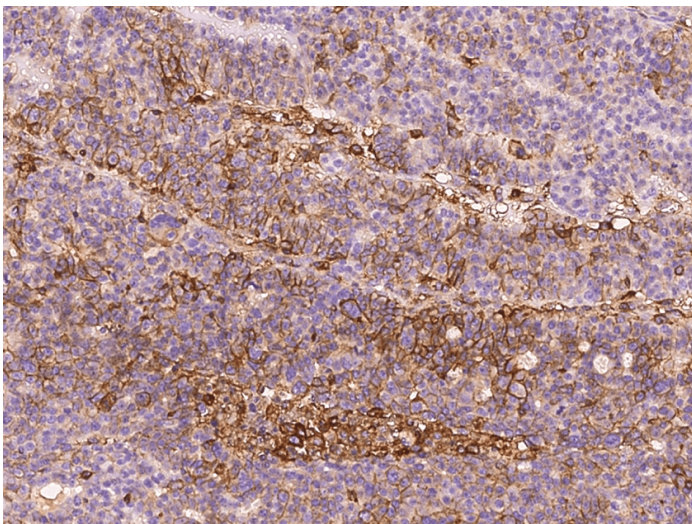
[View Online »](#)



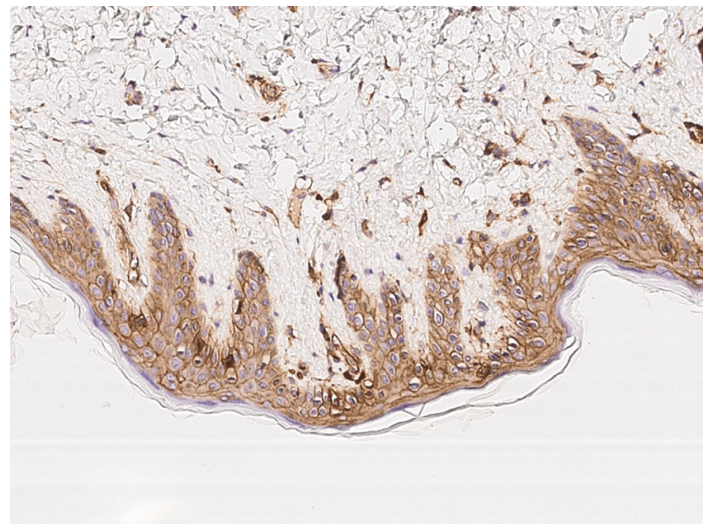
Flow cytometric analysis of Human B2M / Beta-2-microglobulin expression on human whole blood lymphocytes. Cells were stained with purified anti-Human B2M / Beta-2-microglobulin, then a PE-conjugated second step antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of viable lymphocytes.



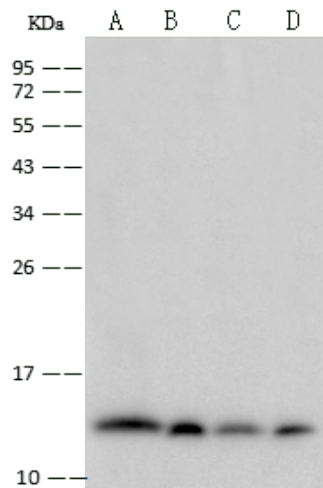
Immunofluorescence staining of B2M in HeLa cells. Cells were fixed with 4% PFA, permeabilized with 0.1% Triton X-100 in PBS, blocked with 10% serum, and incubated with mouse anti-human B2M monoclonal antibody (dilution ratio 1:60) at 4 °C overnight. Then cells were stained with the Alexa Fluor®488-conjugated Goat Anti-mouse IgG secondary antibody (green) and counterstained with DAPI (blue). Positive staining was localized to Cytoplasm and cell membrane.



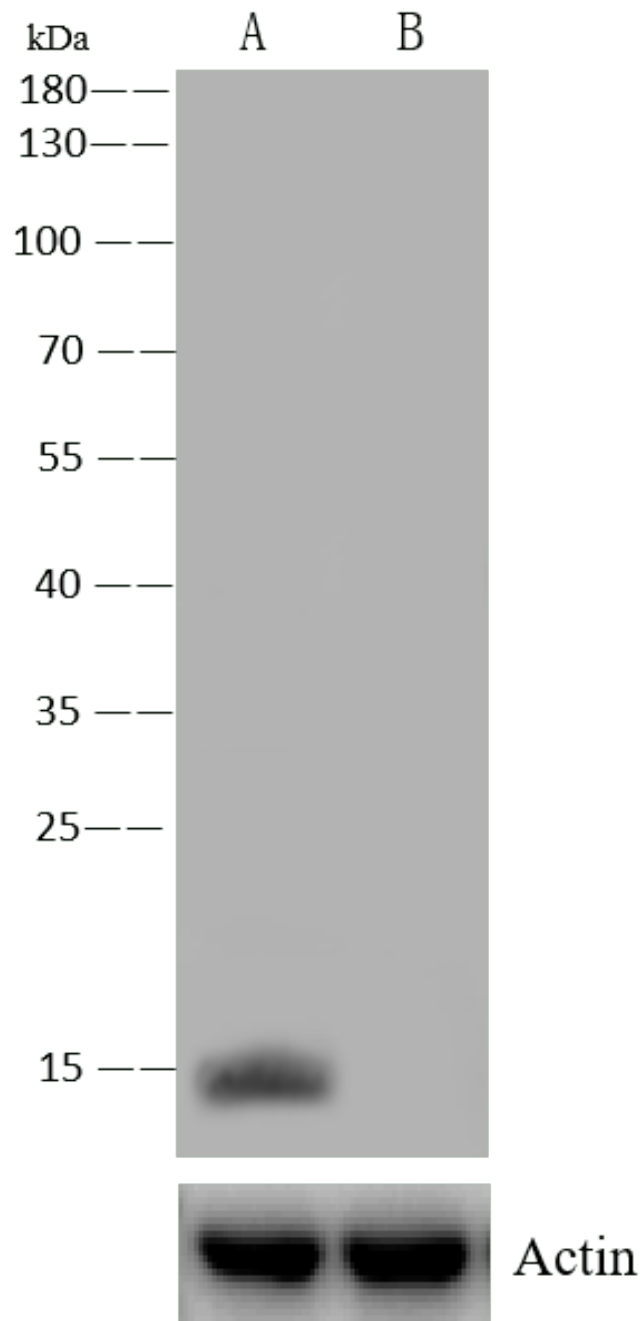
Immunochemical staining of human B2M in human ovarian cancer with mouse monoclonal antibody at 1:60 dilution, formalin-fixed paraffin embedded sections.



Immunochemical staining of human B2M in human skin with mouse monoclonal antibody at 1:60 dilution, formalin-fixed paraffin embedded sections.



Anti-B2M mouse monoclonal antibody at 1:500 dilution. Lane A: A431 Whole Cell Lysate. Lane B: U937 Whole Cell Lysate. Lane C: Raji Whole Cell Lysate. Lane D: HeLa Whole Cell Lysate. Lysates/proteins at 30 µg per lane. Secondary Goat Anti-Mouse IgG (H+L)/HRP at 1/10000 dilution. Developed using the ECL technique. Performed under reducing conditions. Observed band size: 14 kDa



Anti-Beta-2 microglobulin mouse monoclonal antibody at 1:500 dilution. Lane A: HeLa Whole Cell Lysate. Lane B: Beta-2 microglobulin knockout HeLa Whole Cell Lysate. Lysates/proteins at 30 µg per lane. Secondary Goat Anti-Mouse IgG (H+L)/HRP at 1/10000 dilution. Developed using the ECL technique. Performed under reducing conditions. Observed band size: 14 kDa.

Handling & Storage

Handling Protected from light. Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name B2M

Application ELISA, Flow Cytometry, ICC, IF, IHC (PS), WB

Clone 34

Formulation Liquid. 0.2 µm filtered solution in PBS. Preservative free.

Host Mouse

Immunogen Recombinant human Beta-2 microglobulin/B2M protein.

Isotype IgG1

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions Western Blot (1:500-1:1'000)ELISA (1:1'000-1:2'000)Immunohistochemistry (paraffin sections) (1:30-1:100)Immunocytochemistry (1:20-1:100)Immunofluorescence (1:20-1:100)Flow Cytometry (1:25-1:100)ELISA (Detection)(1:1'000-1:10'000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human

Worry-free Guarantee

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



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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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

Last modified: September 17, 2024

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.