

CD63 recombinant monoclonal antibody (001)

This rabbit monoclonal antibody detects CD63 and is used in studies of endosomal trafficking and exosome biology.

CD63 is highly expressed in lysosomes, late endosomes, and extracellular vesicles (EVs). It plays a role in viral infections, cancer metastasis, and immune modulation.

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

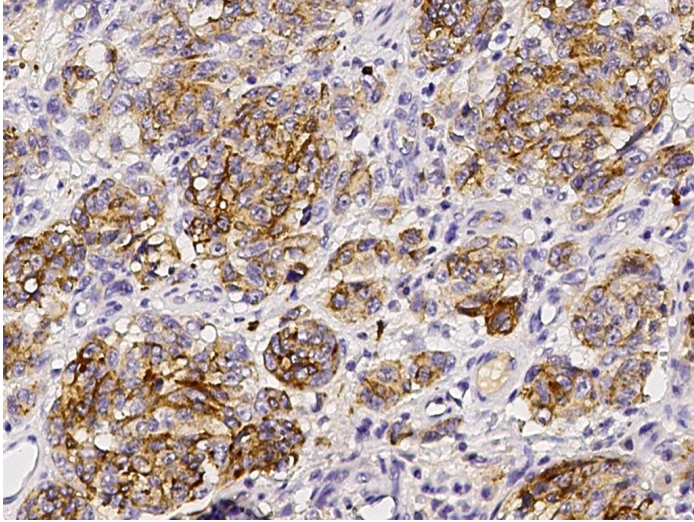
Ordering Information

[Order Online »](#)

ENZ-ABS1068-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunochemical staining of human CD63 in human malignant melanoma with CD63 recombinant monoclonal antibody (ENZ-ABS1068, 1:200, formalin-fixed paraffin embedded sections).

Handling & Storage

Handling Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Tetraspanin-30, Tspan-30

Application IHC (PS)

Clone 001

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

Host Rabbit

Immunogen Recombinant human CD63 / Tspan-30 / Tetraspanin-30 protein.

Isotype IgG

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections) (1:100-1:500)

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-
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)

Last modified: March 28, 2025

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.