

CLIMP-63 monoclonal antibody (G1/296)

The cytoskeleton-linking membrane protein of 62kDa (CLIMP-63) is a marker for the reticular subdomain of the rough endoplasmatic reticulum (ER) and is excluded from the outer nuclear membrane. Concentration of CLIMP-63 into patches may enhance microtubule binding on the cytosolic side and contribute to ER morphology by the formation of a protein scaffold in the lumen of the ER, linking the rough ER to microtubules.

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

Citations: 29

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-ABS669-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	ER protein p63, Cytoskeleton-linking membrane protein of 63kDa
Application	ELISA, Flow Cytometry, ICC, IHC (FS), IP, WB
Clone	G1/296
Formulation	Liquid. In PBS containing 0.1% sodium azide.
Host	Mouse
Immunogen	Human CLIMP-63 (cytoskeleton-linking membrane protein of 63kDa) from Caco-2 cells golgi membrane fraction.
Isotype	IgG2ak
Purity Detail	Protein G-affinity purified.
Source	Purified from hybridoma tissue culture supernatant.
Species Reactivity	Human, Monkey
UniProt ID	Q07065

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: May 29, 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.