

ERp57 polyclonal antibody

ERp57 (also known as Grp58), a member of the Thioredoxin superfamily, is an ER protein with disulfide isomerase activity. It has been shown to increase after oncogenic transformation and to be a component of the MHC Class I peptide-loading complex.

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

Citations: 10

[View Online »](#)

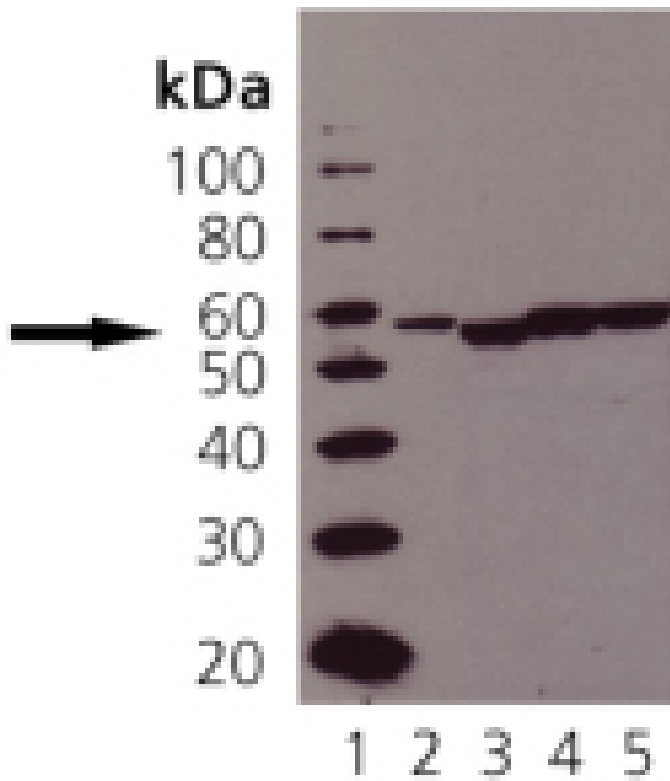
Ordering Information

[Order Online »](#)

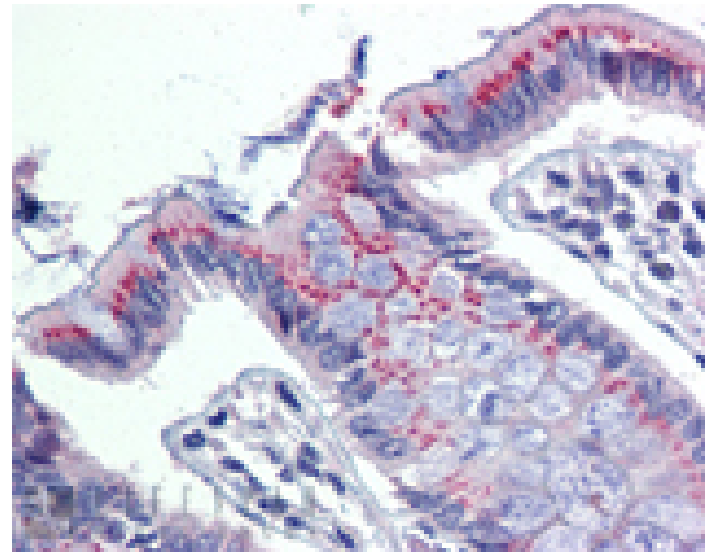
| | |
|---------------|------|
| ADI-SPA-585-D | 50µg |
|---------------|------|

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of ERp57 (Grp58): Lane 1: MW marker, Lane 2: CHO-K1 lysate, Lane 3: Human Liver Microsomes, Lane 4: Rat Liver Microsomes, Lane 5: Mouse Liver Microsomes.



Immunohistochemistry analysis of human colon tissue stained with ERp57, pAb at 10µg/ml.

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 PDIA3, ERp60, ERp61, GRP57, Grp58

Application IHC (PS), WB

Application Notes Detects a band of ~57kDa by Western blot.

Formulation Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

Gene/Protein Identifier NM_017319 (RefSeq)

Host Rabbit

Immunogen Native rat liver ERp57.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:1,000, ECL) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from rabbit serum.

Species Reactivity Bovine, Dog, Guinea pig, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat

UniProt ID P11598

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)