

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 »](#)

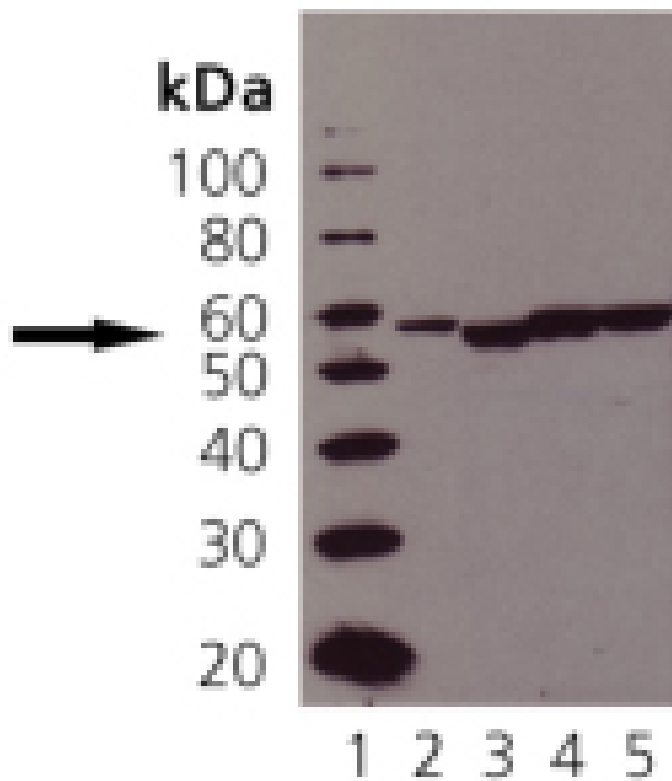
Ordering Information

[Order Online »](#)

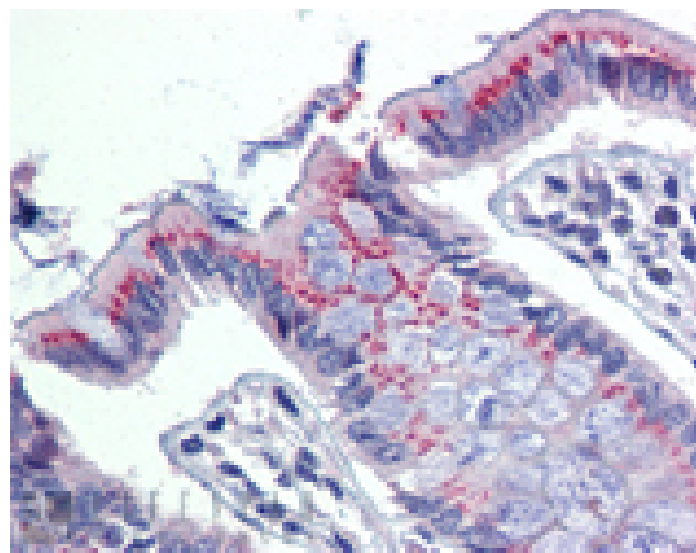
ADI-SPA-585-D	50µg
ADI-SPA-585-F	200µg

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

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



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)