

ERp72 polyclonal antibody

Endoplasmic reticulum protein (ERp) 72, a 72 kDa protein localized to the endoplasmic reticulum, is a member of the Protein Disulfide Isomerase (PDI) family of proteins. Together with other ER resident proteins such as BiP, GRP94, and PDI, they serve as the molecular chaperones for proper folding of newly translocated and glycosylated proteins such as Thyroglobulin (Tg) and Human chorionic gonadotropin (hCG)-beta.

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

Citations: 14

[View Online »](#)

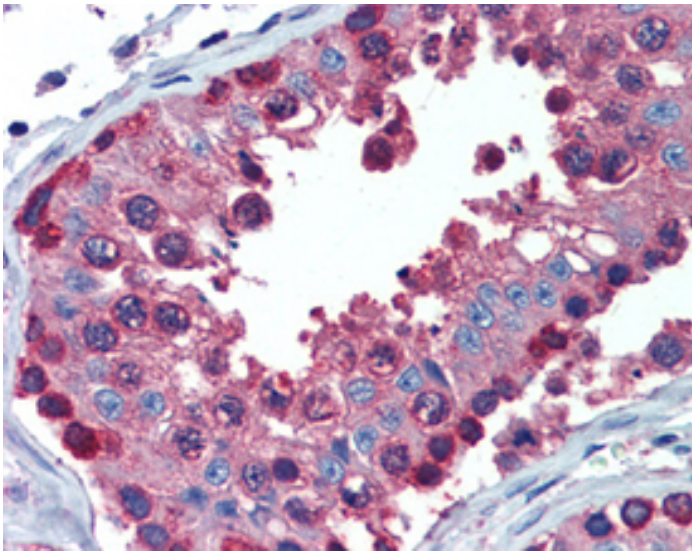
Ordering Information

[Order Online »](#)

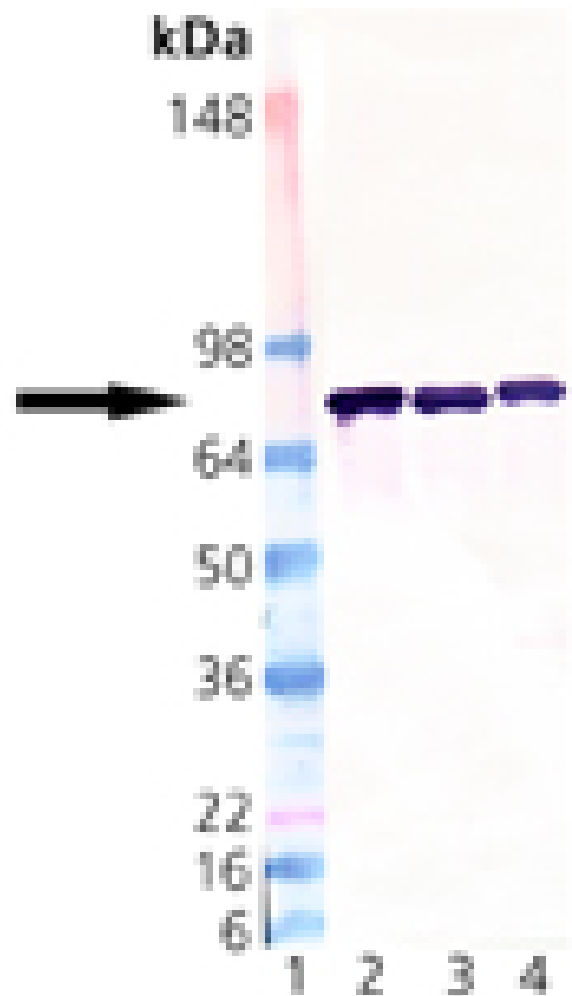
ADI-SPS-720-D	50µg
ADI-SPS-720-F	200µg

Manuals, SDS & CofA

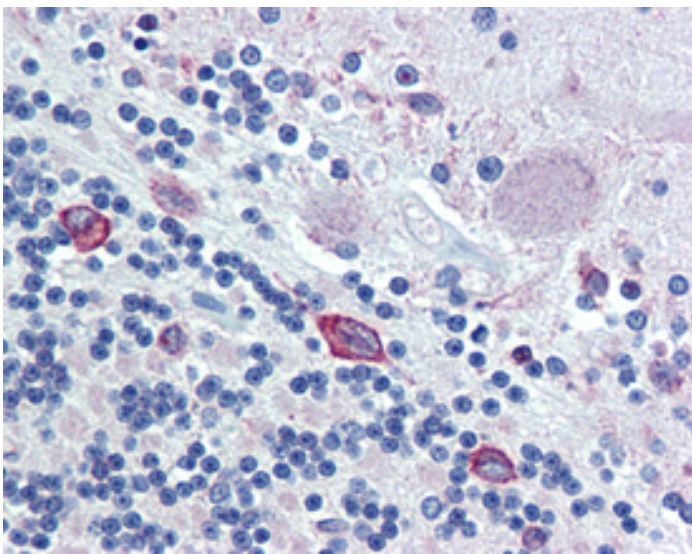
[View Online »](#)



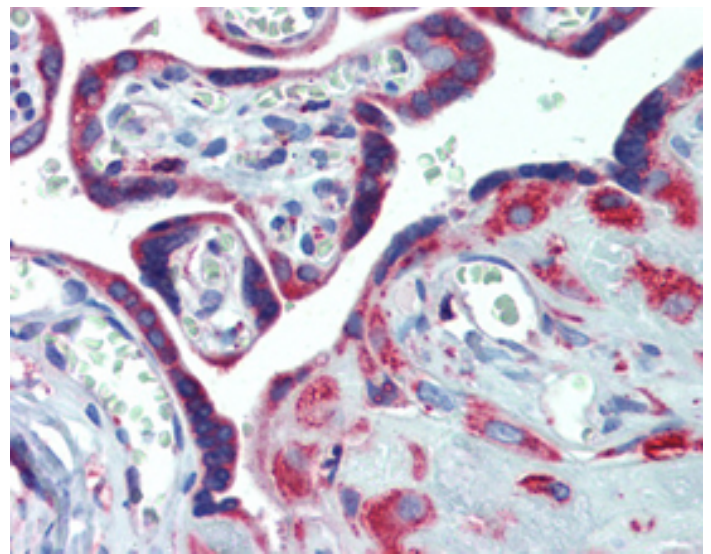
IHC Analysis: Human testis tissue stained with ERp72, pAb at 10µg/ml.



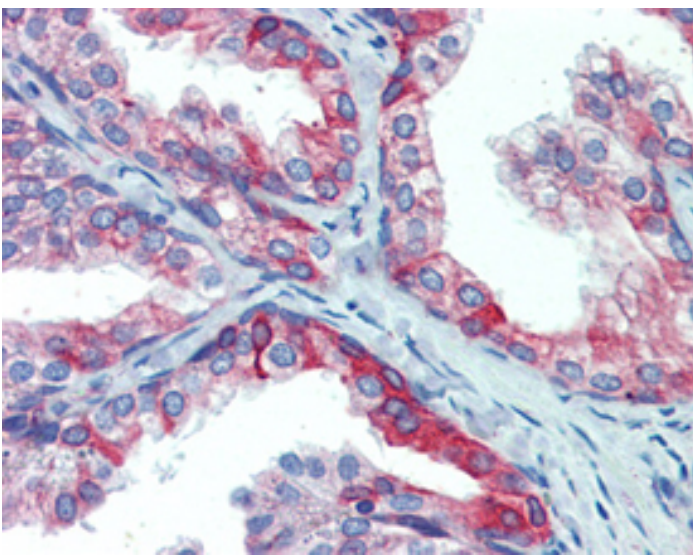
Western blot analysis of ERp72: Lane 1: MW Marker, Lane 2: Rat Liver Microsome Extract, Lane 3: Mouse Liver Microsome Extract, Lane 4: Human Liver Microsome Extract.



IHC Analysis: Human brain, cerebellum tissue stained with ERp72, pAb at 10µg/ml.



IHC Analysis: Human placenta tissue stained with ERp72, pAb at 10µg/ml.



IHC Analysis: Human prostate tissue stained with ERp72, pAb at 10µg/ml.

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

Product Details

Alternative Name	PDIA4, ERP70, protein disulfide isomerase
Application	IF, IHC (PS), WB
Application Notes	Detects a band of ~72kDa by Western blot.
Formulation	Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 623-638 of mouse ERp72.
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:1,000, colorimetric)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, Rabbit, Rat
Technical Info / Product Notes	Recommended by the Human Protein Atlas Organization for IHC (Ensembl No. ENSG00000155660).
UniProt ID	P08003

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)