

Grp94 polyclonal antibody

Grp94 (Glucose-regulated protein 94) is an abundant resident endoplasmic reticulum (ER) luminal stress protein, which together with cytosolic Hsp90 belongs to the Hsp90 family of molecular chaperones. Grp94 expression is upregulated by stress conditions such as glucose starvation and heat shock, which promote protein misfolding or unfolding. In addition to a homeostatic role in protein folding and assembly, Grp94 can function in the intracellular trafficking of peptides from the extracellular space to the MHC class I antigen processing pathway of antigen presentation cells.

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

Citations: 21

[View Online »](#)

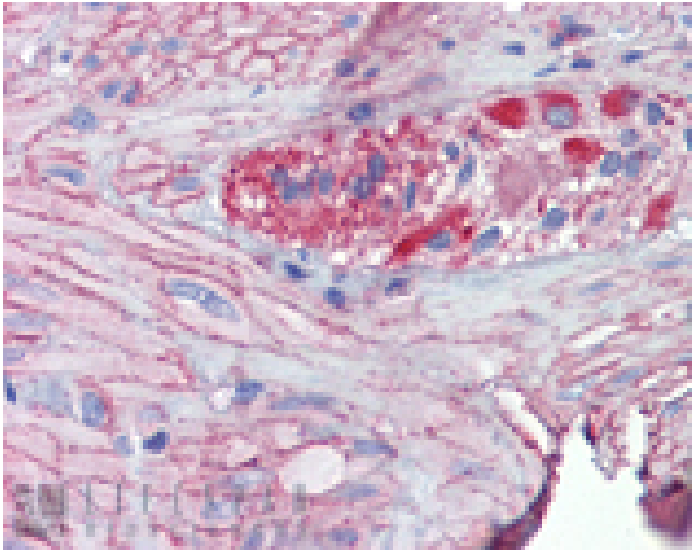
Ordering Information

[Order Online »](#)

ADI-SPA-851-D	50µg
ADI-SPA-851-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of human small intestine tissue stained with Grp94, 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 Endoplasmin, Tra1, Hsp90B1, Gp96

Application ICC, IHC (PS), IP, WB

Application Notes Detects a band of ~94kDa by Western blot.

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

Gene/Protein Identifier NP_035761 (RefSeq)

Host Rabbit

Immunogen Synthetic peptide corresponding to the sequence near the C-terminus of mouse Grp94.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Immunoprecipitation (1:80)Western Blot (1:1,000)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, Human, Mouse, Rat

UniProt ID P08113

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)