

Grp94 (canine), (recombinant)

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.

Citations: 1

[View Online »](#)

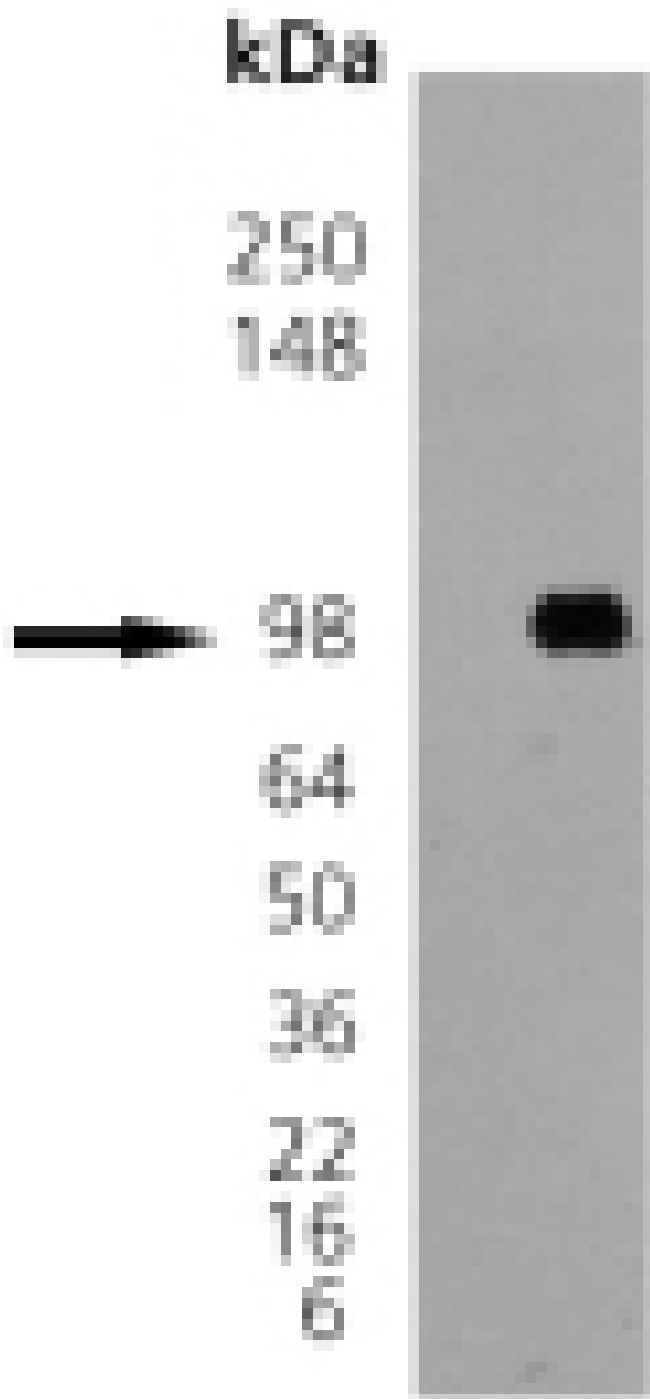
Ordering Information

[Order Online »](#)

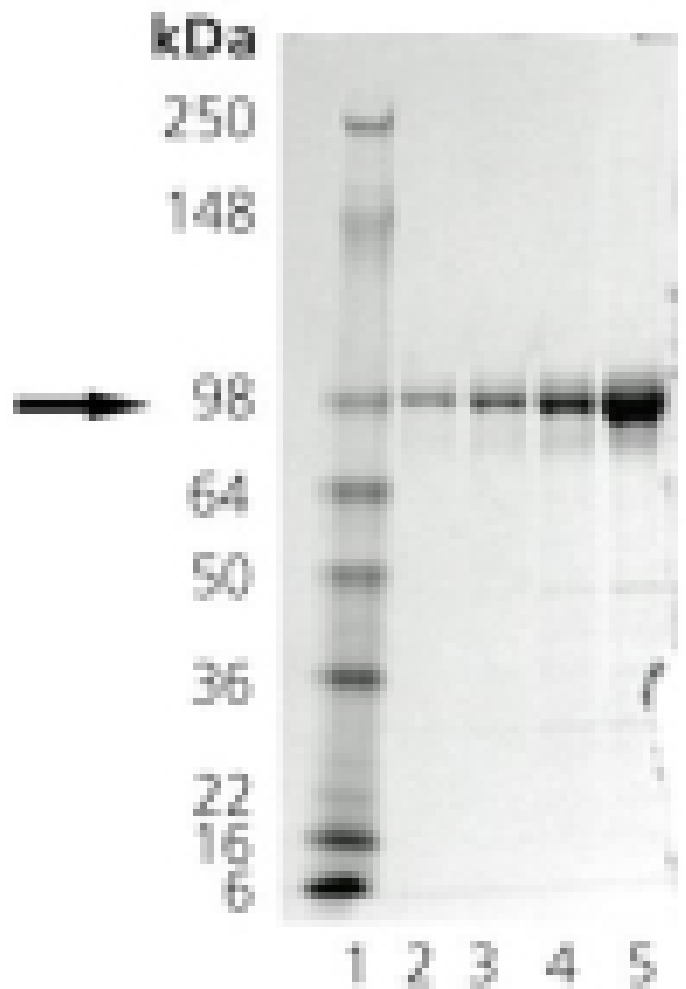
ADI-SPP-766-D	50µg
ADI-SPP-766-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis: 100 ng of Grp94 (canine), (recombinant) probed with Prod. No. ADI-SPA-850.



SDS-PAGE analysis: Lane 1: MW marker, Lane 2: 0.5 µg; Lane 3: 1 µg; Lane 4: 2 µg; Lane 5: 5 µg of Grp94 (canine), (recombinant) detected by Coomassie Stain

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Tra1, Tumor rejection antigen 1, Hsp90B1, Gp96, Heat shock protein 90 kDa beta member 1, Endoplasmin

Application Notes Western blot control.

Formulation Liquid. In 20mM MOPS, pH 7.5, containing 0.5mM EDTA.

MW ~94kDa

Purity ≥90% (SDS-PAGE; Western blot)

Purity Detail Purified by multi-step chromatography.

Source Produced in Sf21 insect cells. Produced in a baculovirus expression system.

UniProt ID P41148



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)