

Grp75/Mortalin (human), (recombinant)

Cells from various organisms increase expression of a class of proteins referred to as heat shock or stress proteins in response to adverse changes in their environment. The members of one class of stress proteins, the Hsp70 family, all bind ATP *in vitro*, but exist within different intracellular compartments. Members include constitutive form Hsc70 within the cytosol/nucleus; inducible form Hsp70 within the cytosol/nucleus/nucleolus; the constitutive glucose-regulated 78 kDa (or BiP) protein within the lumen of the endoplasmic reticulum; and the constitutive glucose-regulated 75 kDa (Grp75) protein within the mitochondrial matrix. Members of the Hsp70 family appear to function as “molecular chaperones, assisting in the folding of other proteins in various intracellular compartments. Grp75 resides in the mitochondrial matrix where it collaborates with Hsp60 in the re-folding of proteins translocated into this organelle. Like its *E. coli* homolog DnaK, Grp75 possesses a cation-dependent ATPase activity considered central to its function as a chaperone.

Citations: 1

[View Online »](#)

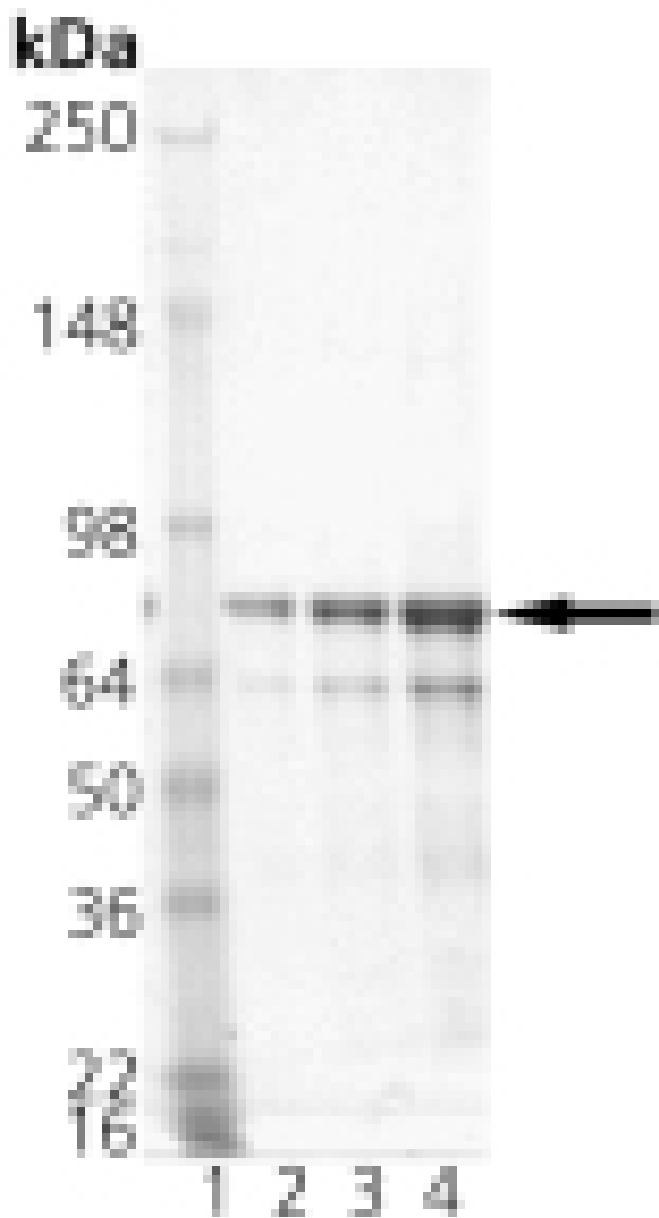
Ordering Information

[Order Online »](#)

ADI-SPP-828-D	50µg
ADI-SPP-828-F	200µg

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis: Lane 1: MW marker, Lane 2: 1 μ g, Lane 3: 2 μ g, Lane 4: 5 μ g of Grp75 (Mortalin) Recombinant Human Protein.



Western Blot analysis: Lane 1: MWM, Lane 2: Grp75 (Mortalin) Recombinant Protein, Lane 3: HeLa Cell Lysate (Heat Shocked) probed with Grp75, pAb.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name MtHsp70, HspA9, Heat shock 70 kDa protein 9

Application Notes Western blot control.

Formulation Liquid. In 50mM TRIS-HCl, pH 7.5, containing 3% glycerol and 150mM sodium chloride.

MW ~75kDa

Purity ≥80% (SDS-PAGE; Western blot)

Purity Detail Purified by multi-step chromatography.

Source Produced in *E. coli*.

UniProt ID P38646



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)