

Grp75/Mortalin monoclonal antibody (30A5)

Grp75 (Glucose regulated 75 kDa protein) is a member of the Hsp70 family. It is localized in the mitochondrial matrix, where, in concert with Hsp60, is thought to participate in the re-folding of proteins translocated into this organelle. Like its *E. coli* homolog DnaK, Grp75 possesses a cation-dependent ATPase activity thought to be central to its function as a chaperone.

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

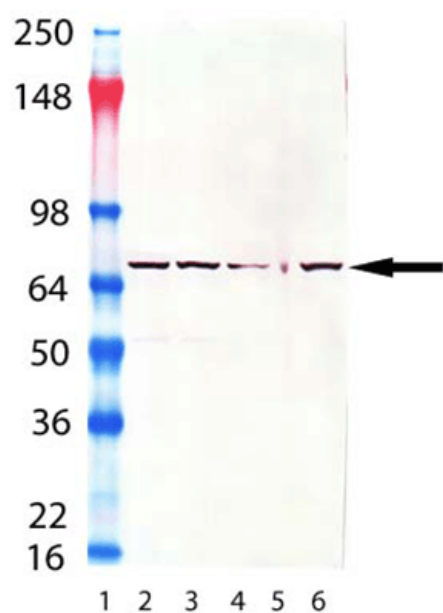
Ordering Information

[Order Online »](#)

ENZ-ABS748-0200	200µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Grp75/Mortalin, mAb (30A5) (Prod. No. ENZ-ABS748): Lane 1: MW marker; Lane 2: HeLa (heat shocked); Lane 3: L-929 (heat shocked); Lane 4: CHO -K1 (heat shocked); Lane 5: GPC-16 (heat shocked); Lane 6: PC-12 (heat shocked).

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	MtHsp70, Hsp70 (mitochondrial), HspA9
Application	IHC, WB
Application Notes	Detects a band of ~75kDa by Western blot.
Clone	30A5
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
GenBank ID	L15189
Host	Mouse
Immunogen	Native human Grp75 (Mortalin).
Isotype	IgG1κ
Purity Detail	Ammonium sulfate precipitation purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (1:50)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 hybridoma tissue culture supernatant.
Species Reactivity	Bovine, Chicken, Dog, Drosophila, Guinea pig, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep, Xenopus
UniProt ID	P38646

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)