

# 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](#).

Citations: 5

[View Online »](#)

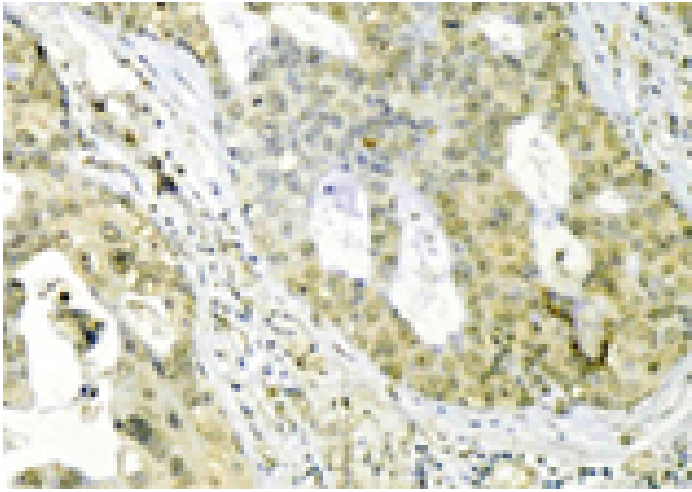
## Ordering Information

[Order Online »](#)

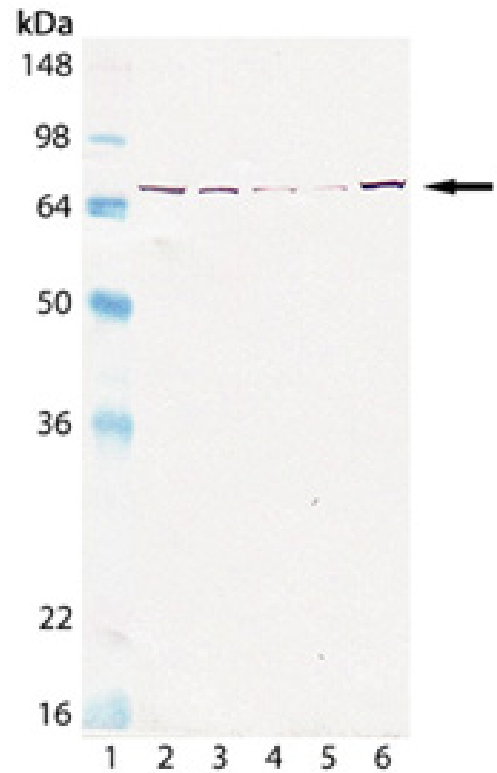
ADI-SPS-825-D	50µl
ADI-SPS-825-F	200µl

## Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of human breast cancer tissue immunohistochemically stained using Grp75 mAb (30A5).



Western blot analysis of Grp75/Mortalin, mAb (30A5) (Prod. no. ADI-SPS-825): 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 IgG1

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 mouse ascites.

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](#)

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)