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

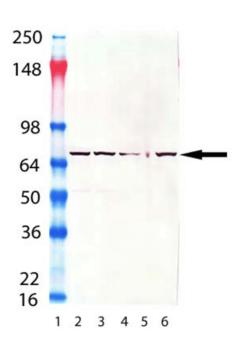
Order Online »

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 NotesDetects 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).

lgG1κ

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

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

ENZO LIFE SCIENCES, European Sales Office ENZO LIFE SCIENCES Phone: 800.942.0430 (ELS) AG Phone: +41 61 926 8989 usa@enzolifesciences.com

eu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 fr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 de@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-uk@enzolifesciences.com