

# HSP70/HSP72 monoclonal antibody (C92F3A-5) (AP conjugate)

The Hsp70 family of heat shock proteins contains multiple homologs ranging in size from 66-78 kDa, and are the eukaryotic equivalents of the bacterial DnaK. The most studied Hsp70 members include the cytosolic stress-induced Hsp70 (Hsp72), the constitutive cytosolic Hsc70 (Hsp73), and the ER-localized BiP (Grp78). Hsp70 family members contain highly conserved N-terminal ATP-ase and C-terminal protein binding domains. Binding of peptide to Hsp70 is assisted by Hsp40, and stimulates the inherent ATPase activity of Hsp70, facilitating ATP hydrolysis and enhanced peptide binding. Hsp70 nucleotide exchange and substrate binding coordinates the folding of newly synthesized proteins, the re-folding of misfolded or denatured proteins, coordinates trafficking of proteins across cellular membranes, inhibits protein aggregation, and targets the degradation of proteins via the proteasomal pathway.

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

Citations: 6

[View Online »](#)

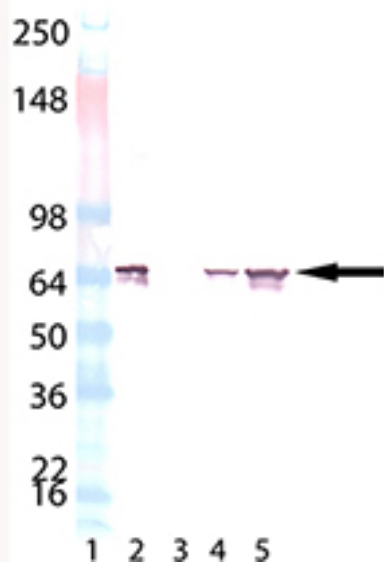
## Ordering Information

[Order Online »](#)

ADI-SPA-810AP-D	50µg
ADI-SPA-810AP-F	200µg

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of HSP70/HSP72, mAb (C92F3A-5) (AP conjugate) (Prod. No. ADI-SPA-810AP): Lane 1: MW marker, Lane 2: HSP70/HSP72 (human), (recombinant) (Prod. No. ADI-NSP-555), Lane 3: HSC70/HSP73 (bovine), (recombinant) (Prod. No. ADI-SPP-751) (Negative Control), Lane 4: HeLa (Prod. No. ADI-LYC-HL100), Lane 5: HeLa (heat shocked) (Prod. No. ADI-LYC-HL101).

# 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                      Hsp70, HspA1A, Heat shock protein 70, HspA1B, Hsp72

Application                      ELISA, WB

Clone                      C92F3A-5

Formulation                      Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

GenBank ID                      M11717

Gene/Protein Identifier                      NP\_005336.3 (RefSeq), NM\_005345 (RefSeq), 3303 (Entrez GeneID), 140550 (OMIM)

Host                      Mouse

Immunogen                      Native human Hsp70.

Isotype                      IgG1

Purity Detail                      Protein G affinity purified.

Recommendation Dilutions/Conditions                      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 ascites.

Species Reactivity                      Bovine, Chicken, Dog, Fish, Human, Mouse, Rat, Teal

UniProt ID

P0DMV8 (HSPA1A), P0DMV9 (HSPA1B)

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)