

# 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 »](#)

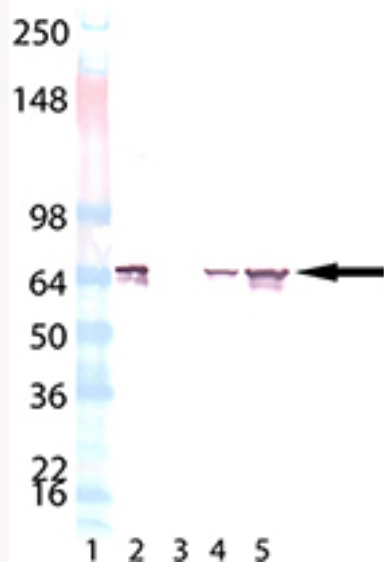
## Ordering Information

[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** Western Blot (1:1,000, colorimetric) Suggested dilutions/conditions may not be available  
**Dilutions/Conditions** 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



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZOLIFE (HSPA1A), RODM9 (HSPA1B)  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
ENZOLIFE (HSPA1A), RODM9 (HSPA1B)  
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)

**Worry-free Guarantee**

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

For local distributors and detailed product information visit us online: [enzo.com](http://enzo.com) All product names, logos, and brands are the property of their respective owners.