

HSC70/HSP73 monoclonal antibody (1B5) (biotin 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 ATPase 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: 7

[View Online »](#)

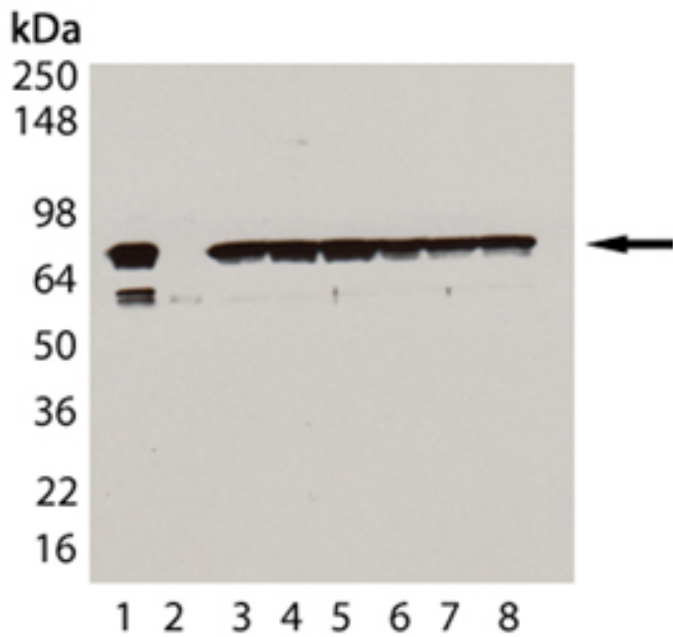
Ordering Information

[Order Online »](#)

ADI-SPA-815B-D	50µg
ADI-SPA-815B-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of HSC70/HSP73, mAb (1B5) (biotin conjugate) (Prod. No. ADI-SPA-815B): Lane 1: HSC70/HSP73 (bovine), (rec) (Prod. No. ADI-SPP-751); Lane 2: HSP70/HSP72 (human), (recombinant) (Prod. No. ADI-NSP-555); Lane 3: 3T3 (heat shocked) (Prod. No. ADI-LYC-3T101); Lane 4: PC-12 (heat shocked) (Prod. No. ADI-LYC-PC101); Lane 5: CHO-K1 (heat shocked); Lane 6: GPC-16 (heat shocked); Lane 7: HeLa (heat shocked) (Prod. No. ADI-LYC-HL101); Lane 8: Vero (heat shocked).

Handling & Storage

Use/Stability	+4°C for short term storage; -20°C for long term storage; Avoid Freeze/Thaw cycles.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Heat shock protein 70, Hsc70, Hsp73
Application	Electron microscopy, ELISA, Flow Cytometry, ICC, IHC, IP, WB
Application Notes	Detects a band of ~73kDa by Western blot.
Clone	1B5
Formulation	Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.
GenBank ID	M34561
Host	Rat
Immunogen	Native hamster Hsc70 (Hsp73).
Isotype	IgG2b
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:1,000, ECL)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, Guinea pig, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep

UniProt ID

P19378

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)