

HSC70/HSP70

monoclonal antibody

(N27F3-4)

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](#).

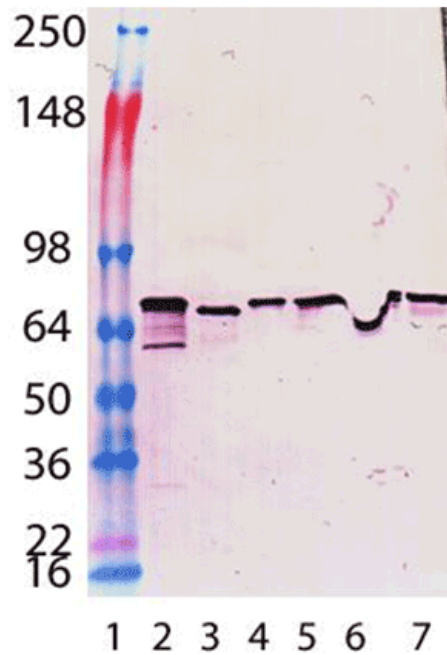
Ordering Information

[Order Online »](#)

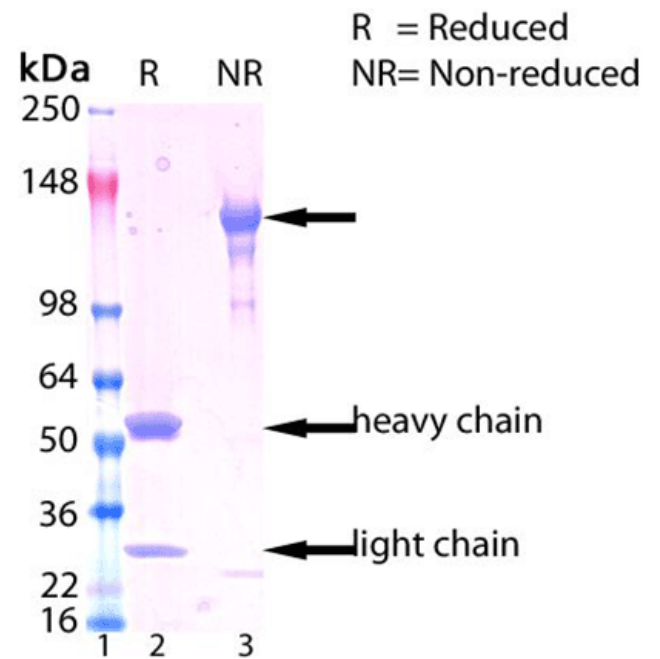
ENZ-ABS716-0050	50µg
ENZ-ABS716-0200	200µg

Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis of ENZ-ABS716: Lane 1: MW Marker, Lane 2: HSC70/HSP73 (bovine), (recombinant) (Prod. No. ADI-SPP-751), Lane 3: HSP70/HSP72 (human), (recombinant) (Prod. No. ADI-NSP-555), Lane 4: 3T3 lysate, Lane 5: 3T3 (heat shocked), Lane 6: HeLa lysate, Lane 7: PC-12 lysate.



SDS-PAGE analysis of ENZ-ABS716: Lane 1: MW Marker, Lane 2 and Lane 3: HSC70/HSP70 monoclonal antibody (N27F3-4), resolved under reduced and non-reduced condition.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Hsc70/Hsp73, Hsp70/Hsp72
Application	Flow Cytometry, IHC (PS), IP, WB
Application Notes	Detects a band of ~72-73kDa by Western blot.
Clone	N27F3-4
Formulation	Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.
GenBank ID	M11717 (HSP70), Y00371 (HSC70)
Host	Mouse
Immunogen	Purified HSC70 (Hsp73) and HSP70 (Hsp72) isolated from human HeLa cells.
Isotype	IgG1κ
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Flow Cytometry (1:100)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	Beluga, Bovine, Chicken, Dog, Fish, Guinea pig, Hamster, Human, Monkey, Mouse, Plant, Porcine, Rabbit, Rat, Sheep, Xenopus
UniProt ID	P0DMV8 (HSP70/HSP72), P11142 (HSC70/HSP73)

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

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

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

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

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