

HSP70 (fish)

polyclonal antibody

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.

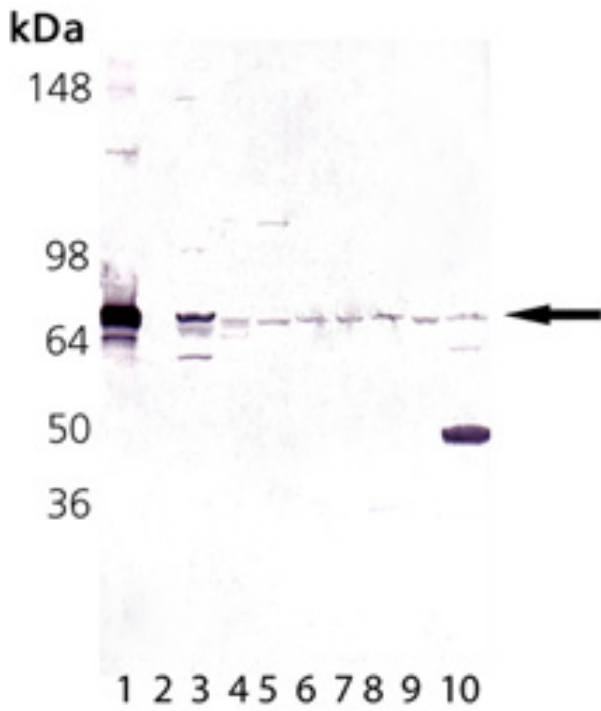
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 2

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Western Blot Analysis of Hsp70 (fish), pAb (Prod. No. ADI-SPA-763): Lane 1: HSP70/HSP72 (Chinook Salmon) (recombinant) (Prod. No. ADI-SPP-763), Lane 2: Dnak (E.Coli) (recombinant) (Prod. No. ADI-SPP-630), Lane 3: RTG-2 Cell Lysate, Lane 4: HeLa Cell Lysate (Prod. No. ADI-LYC-HL100), Lane 5: L-929 Cell Lysate, Lane 6: Rat-2 Cell Lysate, Lane 7: MDBK Cell Lysate, Lane 8: MDCK Cell Lysate, Lane 9: Drosophila Cell Lysate, Lane 10: Chicken Heart Lysate

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 WB

Application Notes Detects a band of ~70kDa by Western blot.

Formulation Liquid. Whole rabbit serum.

Host Rabbit

Immunogen Recombinant salmon Hsp70.

Purity Detail Antiserum.

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 Whole rabbit serum.

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

UniProt ID Q91233

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)

Last modified: May 29, 2024