HSP70 (fish) polyclonal antibody

The Hsp70 family of heat shock protiens 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 »

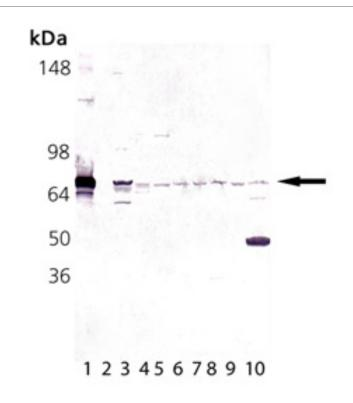
Ordering Information

Order Online »

ADI-SPA-763-D	50µI
ADI-SPA-763-F	200μΙ

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 NotesDetects 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 ReactivityBovine, Chicken, Dog, Drosophila, Fish, Human, Mouse,

Rat

UniProt ID Q91233

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

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