HSC70 (plant) monoclonal antibody (5B7)

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: 9

View Online »

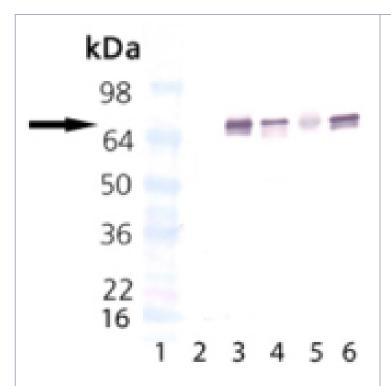
Ordering Information

Order Online »

ADI-SPA-817-D	50µg
ADI-SPA-817-F	200μg

Manuals, SDS & CofA

View Online »



Western blot analysis of HSC70: Lane 1: MW marker, Lane 2: DnaK Recombinant E. coli Protein, Lane 3: Banana Lysate, Lane 4: Orange Lysate, Lane 5: Carrot Lysate, Lane 6: Cytosolic wheat HSC70.

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 Heat shock protein 70, Hsc70, Hsp73

Application IP, WB

Application NotesDetects a band of ~70kDa by Western blot.

Clone 5B7

Formulation Liquid. In PBS containing 50% glycerol and 0.09% sodium

azide.

Host Mouse

Immunogen Recombinant spinach Hsc70 (Hsp73).

lgG1

Purity Detail Protein G affinity purified.

Source Purified from mouse ascites.

Species Reactivity Plant

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

France
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