HSP70/HSP72 monoclonal antibody (N6F3-6.140)

The 70 kDa heat shock protein Hsp70 belongs to the Hsp70 family of highly-related protein isoforms ranging in size from 66 kDa to 78 kDa. Hsc70 shares close biochemical and biological ties to Hsp70, and also belongs to the Hsp70 family. These proteins include cognate members found within major intracellular compartments and highly inducible isoforms predominantly cytoplasmic or nuclear in distribution. Members of the Hsp70 family function as molecular chaperones involved in such cellular functions as protein folding, transport, maturation and degradation, operating in an ATP-dependent manner. The molecular chaperones of the Hsp70 family recognize and bind to nascent polypeptide chains or partially folded intermediates of proteins, preventing their aggregation and misfolding, and the binding of ATP triggers a critical conformational change leading to the release of the bound substrate protein. Data demonstrates that with a ubiquitin-like domain at its amino terminus and its association with the 26S proteosome in HeLa cells, Bag-1 modulates the chaperone activity of Hsc70 and Hsp70. These findings reveal Bag-1's role as a physical link between the Hsc70/Hsp70 chaperone system and the proteasome. Experimental data also shows that the ATPase domain and the substrate binding domain of Hsp70 (or Hsc70) cooperate to form a co-chaperonechaperone complex with the synaptic vesicle cysteine string protein (csp), essential for normal neurotransmitter release.

This antibody is covered by our Worry-Free Guarantee.

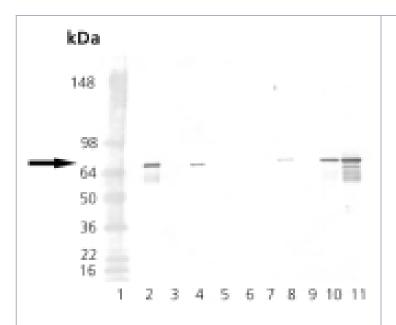
Ordering Information

Order Online »

ADI-SPA-8136-D 50μg

Manuals, SDS & CofA

View Online »



Western blot analysis of HSP70 mAb (N6F3-6.140):
Lane 1: MW marker, Lane 2: HSP70 Recombinant
Human Protein, Lane 3: HSP70-A1 Recombinant
Mouse Protein, Lane 4: DnaK Recombinant E. coli
Protein, Lane 5: HSC70 Recombinant Bovine Protein,
Lane 6: HSC70 (ATPase Fragment) Recombinant
Bovine Protein, Lane 7: HSP70 (HSP72) Recombinant
Rat Protein, Lane 8: HSP70B' Recombinant Human
Protein, Lane 9: HSP70 Recombinant Chinook Salmon
Protein, Lane 10: HeLa, Lane 11: HeLa, Heat Shocked.

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.

Clone N6F3-6.140

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

azide.

GenBank ID M11717

Gene/Protein Identifier NP_005336.3 (RefSeq), NM_005345 (RefSeq), 3303

(Entrez GeneID), 140550 (OMIM)

Host Mouse

Immunogen Native human Hsp70.

lgG1

Purity Detail Protein G affinity purified.

Recommendation Dilutions/ConditionsWestern Blot (1:1,000 colorimetric, 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 hybridoma tissue culture supernatant.

Species Reactivity

UniProt ID

P0DMV8 (HSPA1A), P0DMV9 (HSPA1B)

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

E. coli, Human

Last modified: May 29, 2024



ENZO LIFE SCIENCES,

Phone: 800.942.0430

info-