HSP70-A1 (mouse), (recombinant)

Recombinant mouse HSP70-A1 that functions as a molecular chaperone, assisting protein folding and preventing aggregation.

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.

Citations: 3

View Online »

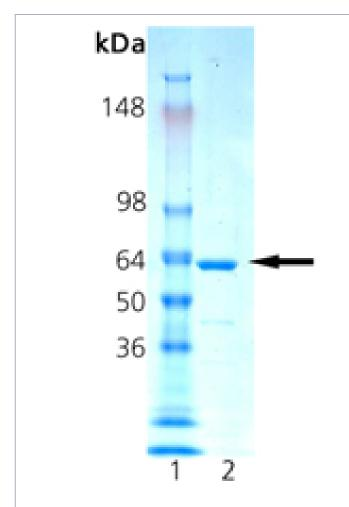
Ordering Information

Order Online »

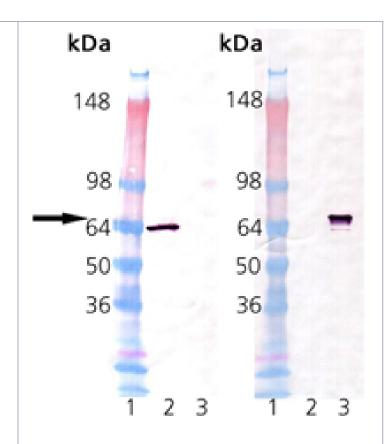
ADI-SPP-502-D	50µg
ADI-SPP-502-F	200μg

Manuals, SDS & CofA

View Online »



SDS-PAGE Analysis: Lane 1: MWM; Lane 2: 1.0 μg of purified Mouse Hsp70-A1 Protein.



Western Blot Analysis: Lane 1: MWM; Lane 2: 100 ng Mouse Hsp70-A1 Protein (Prod. No. ADI-SPP-502); Lane 3: 50 ng E. coli DnaK Protein (Prod. No. ADI-SPP-630); Left: probed with Hsp70 mAb (Prod. No. ADI-SPA-810) at 1.0 μ g/ml. Right: probed with DnaK mAb (Prod. No. ADI-SPA-880) at 1.0 μ g/ml.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Heat shock protein 70

Application NotesATPase activity assay (positive). Western blot control.

Formulation Liquid. In Dulbecco's PBS, pH 7.5, containing 5.0mM

glutathione.

MW ~70kDa (observed)

Purity ≥85% (SDS-PAGE; Western blot)

Purity Detail Purified by multi-step chromatography.

Source Produced in *E. coli*.

UniProt ID Q61696 (Hspa1a), P17879 (Hspa1b)

Last modified: October 9, 2025

