HSP70/HSP72 (Chinook salmon), (recombinant)

Recombinant Chinook salmon HSP70/HSP72, a 70 kDa molecular chaperone that assists in

FINE TO KAP TO PER PROPERTY OF THE PROPERTY OF highly-related protein isoforms ranging in size from 66 kDa to 78 kDa. **prevents aggregation in an ATP-dependent** Hsc70 shares close biochemical and biological ties to Hsp70, and also belongs 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.

Citations: 3

View Online »

Ordering Information

Order Online »

ADI-SPP-763-F

200µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Heat shock protein 70, HspA1A, HspA1B

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

Formulation Liquid. In Dulbecco's PBS.

MW ~70kDa

Purity ≥90% (SDS-PAGE; Western blot)

Purity Detail Purified by multi-step chromatography.

Source Produced in *E. coli*.

Last modified: October 9, 2025



eu@enzolifesciences.com