

HSP70 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # H34-54H

Lot # R087-1

Product Description

Recombinant full-length human HSP70 was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag. The gene accession number is [NM_005345](#).

Gene Aliases

HSPA1A, HSP72, HSPA1, HSPA1B, HSP70-1

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

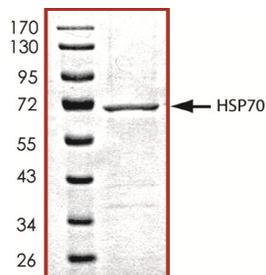
Scientific Background

HSP70 is a member of the heat shock protein family and is synthesized by cells of many organisms in response to stress (1). HSP70 is found mostly but not exclusively in the nucleus of unstressed cells. For several hours after a short heat shock, it is strongly concentrated in nucleoli (2). Nucleoli are transiently damaged by such a heat shock. Their morphology changes and assembly and export of ribosomes are blocked for several hours. HSP70 helps to stabilize the morphological changes of the nucleoli.

References

1. Moran, L A. et al: The major heat-shock protein (hsp70) gene family: related sequences in mouse, Drosophila, and yeast. *Can J Biochem Cell Biol.* 1983 Jun;61(6):488-99.
2. Pelham, H R: Hsp70 accelerates the recovery of nucleolar morphology after heat shock. *EMBO J.* 1984 Dec 20;3(13):3095-100.

Purity



The purity of HSP70 protein was determined to be **>90%** by densitometry. Approx. MW **70kDa**.

Upstream Active Kinases

AKT1	Cat # A16-10G
AKT2	Cat # A17-10G
AKT3	Cat # A18-10G

HSP70 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	H34-54H
Specific Lot Number	R087-1
Purity	>90%
Concentration	0.2 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.