

## HSP40 Protein

Full-length recombinant protein expressed in E. coli cells

**Catalog # H32-54G**

Lot # K208-1

### Product Description

Recombinant full-length human HSP40 was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM\\_006145](#).

### Gene Aliases

DNAJB1; HSPF1

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 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

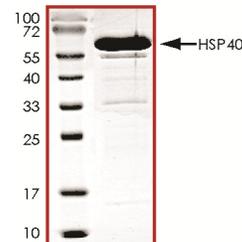
HSP40 is a member of heat shock proteins (HSPs) which are synthesized in cells in response to heat shock and other metabolic stresses and provide a transient state of thermotolerance. HSP40 is localized faintly throughout the cell in non-heat-shocked cells and accumulates in the nuclei and nucleoli in heat-shocked cells (1). The intracellular localization of HSP40 is very similar to that of HSP70. Overexpression of HSP40 reduces aggregate formation and suppresses apoptosis in a neuronal cell model of spinal and bulbar muscular atrophy (inherited neurodegenerative diseases caused by polyglutamine expansion) (2).

### References

- Hattori, H. et al: Intracellular localization and partial amino acid sequence of a stress-inducible 40-kDa protein in HeLa cells. *Cell Struct. Funct.* 1992;17(1):77-86.
- Kobayashi, Y. et al: Chaperones Hsp70 and Hsp40 suppress aggregate formation and apoptosis in cultured neuronal cells expressing truncated androgen receptor protein with expanded polyglutamine tract. *J Biol Chem.* 2000 24;275:8772-8

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

### Purity



The purity of HSP40 was determined to be **>90%** by densitometry. Approx. MW **65kDa**.

### Upstream Active Kinases

MAPKAPK2	Cat # M40-11
MAPKAPK3	Cat # M41-10

## HSP40 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog Number	H32-54G
Specific Lot Number	K208-1
Purity	>90%
Concentration	0.2 µg/µl
Stability	1 yr 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.