

Catalogue #	Aliquot Size
H36-50H-50	50 µg
H36-50H-200	200 µg

HSP90 α Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # H36-50H

Lot # R279-1

Product Description

Recombinant full length human HSP90 α was expressed by baculovirus in Sf9 insect cells using a C-terminal His tag. The gene accession number is [NM_005348](#).

Gene Aliases

HSP90AA1, HSPN, LAP2, HSP86, HSPC1, HSPCA, Hsp89, Hsp90, HSP90N, HSPCAL1, HSPCAL4, FLJ31884

Formulation

Recombinant protein stored in 50mM MOPS, 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.

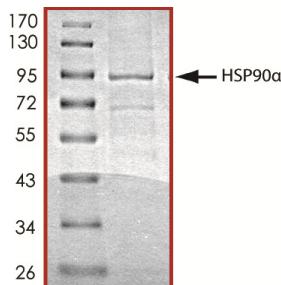
Scientific Background

HSP90 α is a member of the HSP90 family of proteins which are important molecular chaperones involved in signal transduction, cell cycle control, stress management, folding, degradation, and transport of proteins (1). HSP90 is a molecular chaperone that plays a key role in the conformational maturation of oncogenic signaling proteins, including HER2/ERBB2, AKT, RAF1, BCR-ABL and mutated p53. HSP90 inhibitors bind to HSP90, and induce the proteasomal degradation of HSP90 client proteins. HSP90 α is an important mediator of cancer cell invasion and is expressed extracellularly on fibrosarcoma and breast cancer cells where it interacts with MMP2 (2).

References

1. Csermely, P. et al: The 90-kDa molecular chaperone family: structure, function, and clinical applications. *Pharmacol Ther.* 1998 Aug;79(2):129-68.
2. Eustace, B. K. et al: Functional proteomic screens reveal an essential extracellular role for hsp90-alpha in cancer cell invasiveness. *Nature Cell Biol.* 6: 507-514, 2004.

Purity



The purity of HSP90 α was determined to be >85% by densitometry.
Approx. MW 94kDa.

HSP90 α Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	H36-50H
Specific Lot Number	R279-1
Purity	>85%
Concentration	0.2µg/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.