

HSP90 β Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # H36-54H

Lot #G324-1

Product Description

Recombinant full length human HSP90 β was expressed by baculovirus in Sf9 insect cells using a C-terminal His tag. The gene accession number is [NM_007355](#).

Gene Aliases

HSP90AB1, HSPC2, HSPCB, D6S182, HSP90B

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.2mM 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

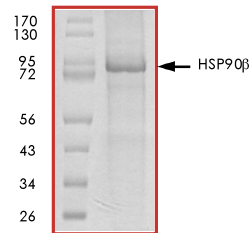
HSP90 β is a member of the HSP90 family of proteins which are important molecular chaperones involved in signal transduction, cell cycle control, stress management, folding, degradation, and transport of proteins (1). HSP90 proteins have been found in a variety of organisms suggesting that they are ancient and conserved. HSP90 binds to client proteins (such as steroid receptors, AKT, Bcr-Abl, Apaf-1, survivin, cyclin dependent kinases) and acts as a molecular chaperone. Failure of HSP90 chaperone activity leads to misfolding of client proteins, which leads to ubiquitination and proteasome degradation, and thus deregulation of cellular homeostasis (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. Georgakis, G.V. et al: Heat-shock protein 90 inhibitors in cancer therapy: 17AAG and beyond. *Future Oncol.* 2005 Apr;1(2):273-81.

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

Purity



The purity was determined to be **>90%** by densitometry.
Approx. MW **91kDa**.

Upstream Active Kinases

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

HSP90 β Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	H36-54H
Specific Lot Number	G324-1
Purity	>90%
Concentration	0.2 $\mu\text{g}/\mu\text{l}$
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.