

Anti-HSP90β

Rabbit Polyclonal Antibody

Catalog # H36-63R

Lot # H200-4

Cited Applications

- Western blot (1:1000)

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the HSP90β protein

Cross Reactivity

- Western blot of human and mouse HSP90β

HSP90β from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to amino acids 252-256 of human HSP90β

Formulation

PBS pH 7.4, 0.02% sodium azide, 50% glycerol

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

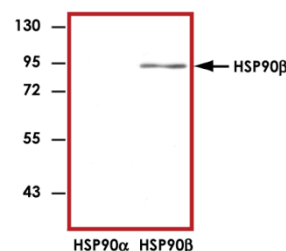
HSP90β is a member of the HSP90 family of proteins which are important molecular chaperones involved in signal transduction, cell cycle control, stress management, and folding, degradation, and transport of proteins (1). HSP90 proteins have been found in a variety of organisms suggesting that they are 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

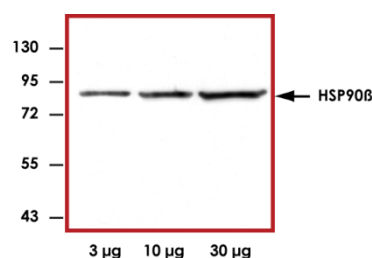
1. Csermely, P. et al: The 90-kDa molecular chaperone family: structure, function, and clinical applications. *Pharmacol Ther.* 1998 Aug;79(2):129-68.
2. Georgakias, G.V. et al: Heat-shock protein 90 inhibitors in cancer therapy: 17AAG and beyond. *Future Oncol.* 2005 Apr;1(2):273-81.

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

Sample Data



Representative western blot with Anti-HSP90β (1:1000) using 20ng of human recombinant HSP90α and HSP90β.



Representative western blot with Anti-HSP90β (1:1000) using 3 µg, 10 µg, and 30 µg of HeLa cell lysate.

Anti-HSP90β

Rabbit Polyclonal Antibody

Catalog Number

H36-63R

Specific Lot Number

H200-4

Purification

Affinity chromatography

Concentration

1.0 µg/µL

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.