HSP90β polyclonal antibody

The 90 kDa molecular chaperone family includes 90 kDa heat shock protein Hsp90 and 94 kDa glucose-regulated protein grp94, both major molecular chaperones of the cytosol and the endoplasmic reticulum. Mammalian cells express inducible Hsp90 alpha and constitutive Hsp90 beta isoforms that are encoded by separate genes. The amino acid sequences of human and yeast Hsp90 alpha are 85% and 90% homologous to those of Hsp90 beta, respectively. All known members of the Hsp90 protein family are highly conserved, especially in the N-terminal and C-terminal regions containing independent chaperone sites with different substrate specificity. These ubiquitous and highly conserved proteins account for 1-2% of all cellular protein in most cells. Hsp90 functions as part of the cell's powerful network of chaperones to fight the deleterious consequences of protein unfolding caused by non-physiological conditions. In the absence of stress, however, Hsp90 provides a necessary component of such fundamental cellular processes as hormone signaling and cell cycle control by serving as a chaperone for many key signaling molecules including steroid receptors, cell cycle kinases involved in signal transduction, and p53. As many of these client proteins are known oncogenes, Hsp90 inhibitors such as 17-AAG, a geldanamycin analog, have been of benefit in the treatment of many cancers.

This antibody is covered by our Worry-Free Guarantee.

Citations: 7

View Online »

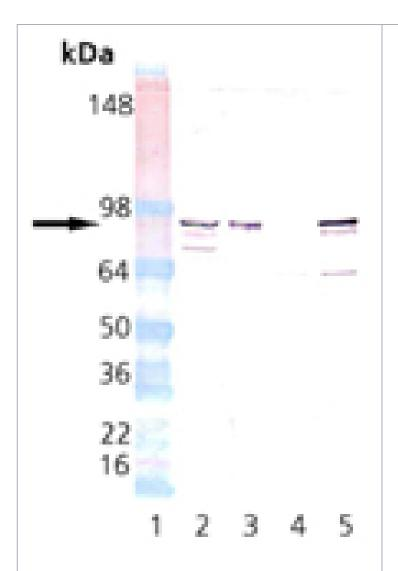
Ordering Information

Order Online »

ADI-SPA-844-050	50µg
ADI-SPA-844-200	200μg

Manuals, SDS & CofA

View Online »



Western blot analysis of HSP90 beta pAb: Lane 1: MW marker, Lane 2: HeLa Cell Lysate, Heat Shocked (Prod. No. ADI-LYC-HL101), Lane 3: HSP90 Native Human Protein (Prod. No. ADI-SPP-770), Lane 4: HSP90 alpha Recombinant Human Protein (Prod. No. ADI-SPP-776), Lane 5: HSP90 beta Recombinant Human Protein (Prod. No. ADI-SPP-777).

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name HSP86, Heat shock protein 90

Application WB

Application NotesDetects a band of ~90kDa by Western blot.

Formulation Liquid. In PBS containing 50% glycerol and 0.09% sodium

azide.

Host Rabbit

Immunogen Recombinant human HSP90β.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/ConditionsWestern Blot (1:1,000, colorimetric)Suggested

dilutions/conditions may not be available for all applications. Optimal conditions must be determined

individually for each application.

Source Purified from rabbit serum.

Species Reactivity Bovine, Dog, Guinea pig, Hamster, Human, Mouse,

Porcine, Rabbit, Sheep

UniProt ID P08238

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com