HSP90α monoclonal antibody (K41009)

The Hsp90 family of heat shock proteins represents one of the most abundantly expressed and highly conserved families of cellular chaperones whose expression can be upregulated under conditions of cellular stress, and includes cytoplasmic (Hsp90 α / β), ER (grp94), and mitochondrial (TRAP1) localized members. Structurally, Hsp90 is characterized by an N-terminal ATP-binding domain, a medial substrate-binding domain, and a C-terminal dimerization motif. Hsp90 dimers function in cooperation with cochaperones (e.g. Hsp40, Hsp70, Hop, p23) to stabilize a multitude of client protein substrates, including steroid hormone receptors, protein kinases, and transcription factors. The essential binding and hydrolysis of ATP by Hsp90 is inhibited by ansamycin drugs (e.g. geldanamycin, 17-AAG) which occupy the N-terminal Hsp90 nucleotide-binding pocket. Many Hsp90 client proteins such as erbB2/Her-2, c-raf, bcr-abl, p53, and hTERT, are members of well characterized oncogenic pathways, making Hsp90 inhibitors useful anticancer agents.

This antibody is covered by our Worry-Free Guarantee.

Citations: 3

View Online »

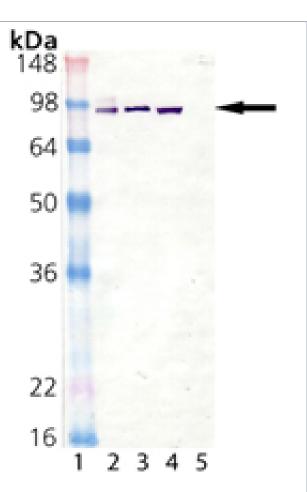
Ordering Information

Order Online »

| ADI-SPA-839-050 | 50µg |
|-----------------|-------|
| ADI-SPA-839-200 | 200μg |

Manuals, SDS & CofA

View Online »



Western blot analysis of HSP90, mAb (K41009): Lane 1: Marker; Lane 2: HSP90 (human), (native) (Prod. No. ADI-SPP-770); Lane 3: HSP90 α (human), (recombinant) (Prod. No. ADI-SPP-776); Lane 4: HeLa (heat shocked, (cell lysate) (Prod. No. ADI-LYC-HL101); Lane 5: HSP90 β (human), (recombinant) (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 90α

Application WB

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

Clone K41009

Formulation Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.

Host Mouse

Immunogen Recombinant human Hsp90α.

Isotype IgG2a

Purity Detail Protein G-affinity purified.

Recommendation Dilutions/Conditions Western Blot (1µg/ml, colorimetric)Suggested

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

individually for each application.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human, Mouse, Rabbit, Rat

UniProt ID P07900

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

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com