

# HSP90 $\alpha$ 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 $\alpha/\beta$ ), 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 »](#)

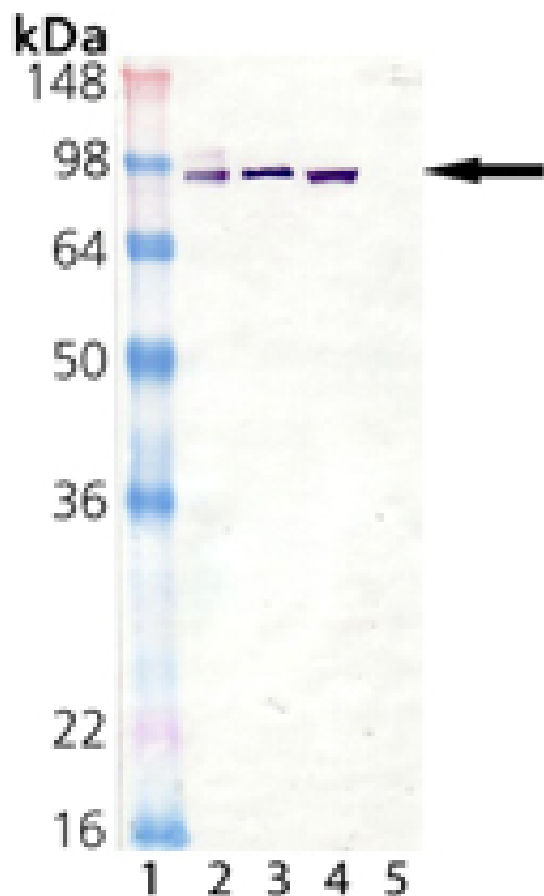
## Ordering Information

[Order Online »](#)

ADI-SPA-839-050	50 $\mu$ g
ADI-SPA-839-200	200 $\mu$ g

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 $\alpha$  (human), (recombinant) (Prod. No. ADI-SPP-776); Lane 4: HeLa (heat shocked, (cell lysate) (Prod. No. ADI-LYC-HL101); Lane 5: HSP90 $\beta$  (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 Notes** Detects 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** Western Blot (1µg/ml, colorimetric) Suggested dilutions/conditions may not be available  
**Dilutions/Conditions** 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



ENZO LIFE SCIENCES,  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 1 42 44 55 55  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

This antibody is covered by our **Worry Free Guarantee**