

HSP90 β monoclonal antibody (K3705)

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: 8

[View Online »](#)

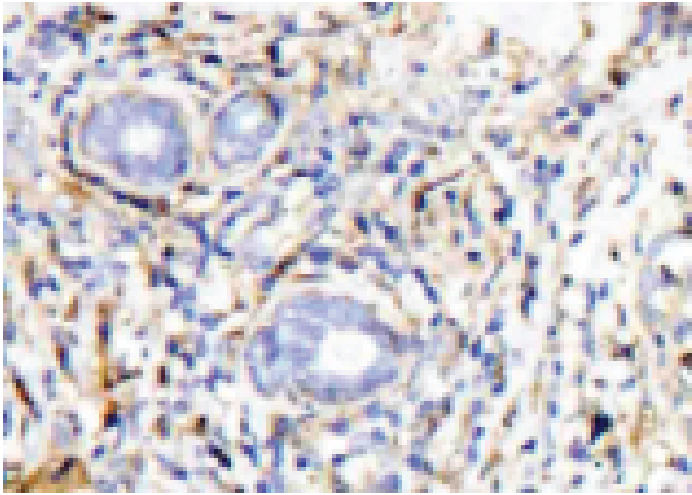
Ordering Information

[Order Online »](#)

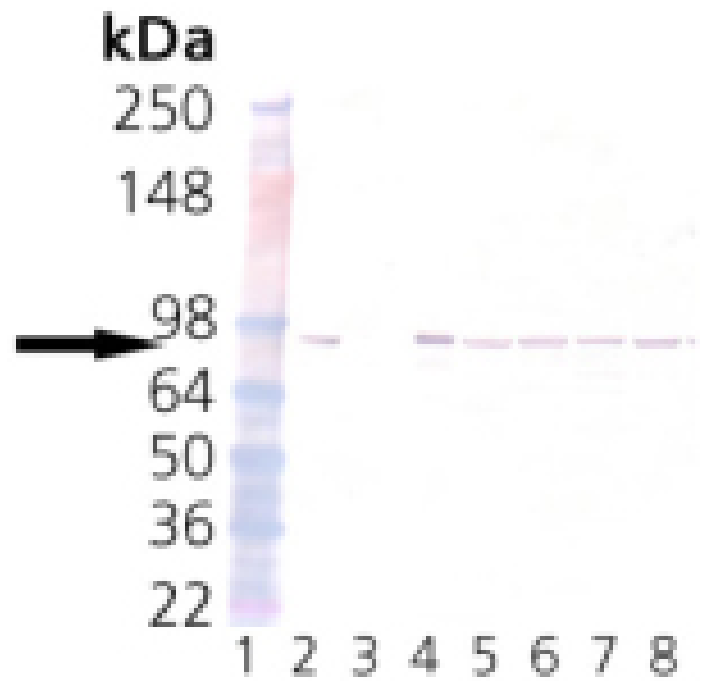
ADI-SPA-842-F	200 μ g
---------------	-------------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of human breast cancer tissue using Hsp90 β mAb (K3705).



Western blot analysis of HSP90 β : Lane 1: MW marker, Lane 2: HSP90 Native Human Protein, Lane 3: HSP90 α Recombinant Human Protein, Lane 4: HSP90 β Recombinant Human Protein, Lane 5: HeLa, Lane 6: PC-12, Lane 7: CHO-K1, Lane 8: 3T3.

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 90β

Application IHC, WB

Application Notes Detects a band of ~90kDa by Western blot.

Clone K3705

Formulation Liquid. In saline, pH 6.9, containing 0.1% sodium azide.

Host Mouse

Immunogen Recombinant human HSP90β.

Isotype IgM

Purity Detail Ammonium sulfate precipitation purified.

Recommendation Dilutions/Conditions Immunohistochemistry (1:50)Western 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 hybridoma tissue culture supernatant.

Species Reactivity Bovine, Chicken, Dog, Guinea pig, Hamster, Human, Mouse, Porcine, Rat, Sheep

UniProt ID P08238



ENZO LIFE SCIENCES,
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 1 42 44 55 55
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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