

HSP90 β monoclonal antibody (K3701)

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: 4

[View Online »](#)

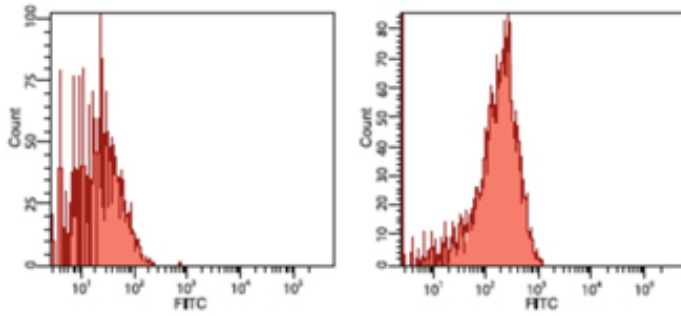
Ordering Information

[Order Online »](#)

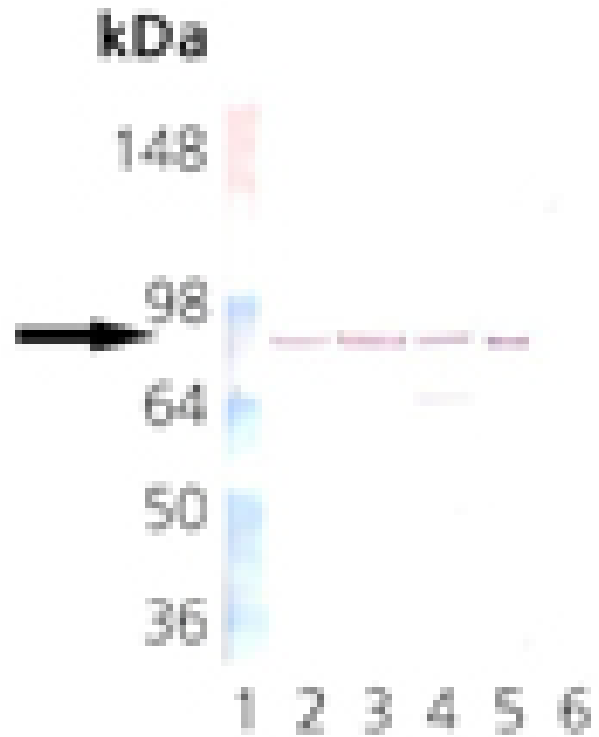
ADI-SPA-843-D	50 μ g
ADI-SPA-843-F	200 μ g

Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis human colon cancer Caco-2 cells analyzed by flow cytometry using isotype control antibody (left) or Hsp90 β mAb (K3701) (right).



Western blot analysis: Lane 1: MW marker, Lane 2: HeLa, Lane 3: PC-12, Lane 4: CHO-K1, Lane 5: HSP90 Native Protein (Prod. No. ADI-SPP-770), Lane 6: HSP90 alpha Recombinant Protein (Prod. No. ADI-SPP-776) (negative control) , probed with HSP90 β mAb (K3701)

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	HSP84, Heat shock protein 90β
Application	Flow Cytometry, WB
Application Notes	Detects a band of ~90kDa by Western blot.
Clone	K3701
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	Flow Cytometry (10µg/ml)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, Dog, Guinea pig, Hamster, Human, Mouse, Porcine, Rabbit, Rat, Sheep
UniProt ID	P08238

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



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

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

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

France
Phone: +33 472 440 655
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