HSP90β 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, bcrabl, 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 »

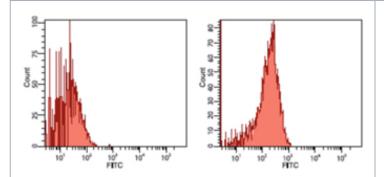
Ordering Information

Order Online »

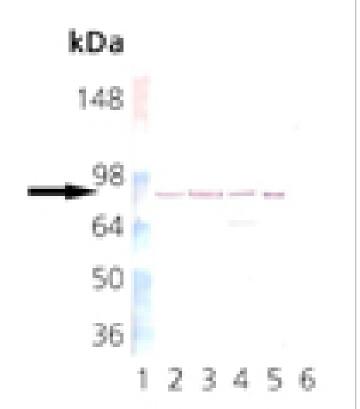
ADI-SPA-843-D	50µg
ADI-SPA-843-F	200μg

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: HeL a, 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 NotesDetects 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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com