

# HSP90 $\beta$ polyclonal antibody (PE conjugate)

The 90 kDa molecular chaperone family includes 90 kDa heat shock protein Hsp90 and 94 kDa glucose-regulated protein grp94, both major molecular chaperones of the cytosol and the endoplasmic reticulum. Mammalian cells express inducible Hsp90 $\alpha$  and constitutive Hsp90 $\beta$  isoforms that are encoded by separate genes. The amino acid sequences of human and yeast Hsp90 $\alpha$  are 85% and 90% homologous to those of Hsp90 $\beta$ , respectively. All known members of the Hsp90 protein family are highly conserved, especially in the N-terminal and C-terminal regions containing independent chaperone sites with different substrate specificity. These ubiquitous and highly conserved proteins account for 1-2% of all cellular protein in most cells. Hsp90 functions as part of the cell's powerful network of chaperones to fight the deleterious consequences of protein unfolding caused by non-physiological conditions. In the absence of stress, however, Hsp90 provides a necessary component of such fundamental cellular processes as hormone signaling and cell cycle control by serving as a chaperone for many key signaling molecules including steroid receptors, cell cycle kinases involved in signal transduction, and p53. As many of these client proteins are known oncogenes, Hsp90 inhibitors such as 17-AAG, a geldanamycin analog, have been of benefit in the treatment of many cancers.

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

Citations: 5

[View Online »](#)

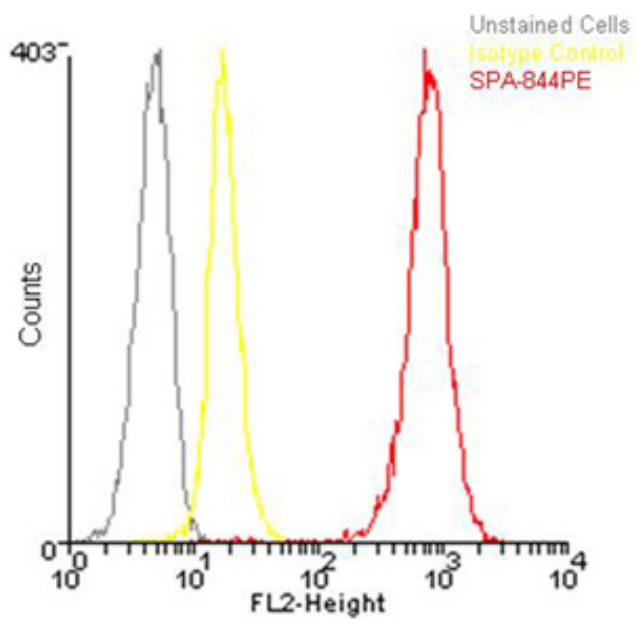
## Ordering Information

[Order Online »](#)

ADI-SPA-844PE-050	50 $\mu$ g
ADI-SPA-844PE-200	200 $\mu$ g

## Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of  $10^6$  Jurkat cells using Hsp90 $\beta$ , pAb (PE conjugate) at a concentration of 50 $\mu$ g/ml.

## Handling & Storage

**Handling** Avoid freeze/thaw cycles. Protect from light.

**Long Term Storage** +4°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** HSP84, Heat shock protein 90β

**Application** Flow Cytometry

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

**Formulation** Liquid. In PBS containing 0.09% sodium azide.

**Host** Rabbit

**Immunogen** Recombinant human HSP90β.

**Purity Detail** Protein A affinity purified.

**Recommendation Dilutions/Conditions** Flow Cytometry (1:100) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

**Source** Purified from rabbit serum.

**Species Reactivity** Bovine, Dog, Guinea pig, Hamster, Human, Mouse, Porcine, Rabbit, Sheep

**UniProt ID** P08238

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



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

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

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

France  
Phone: +33 472 440 655  
[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)