

HSP90 monoclonal antibody (AC88) (PE conjugate)

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 »](#)

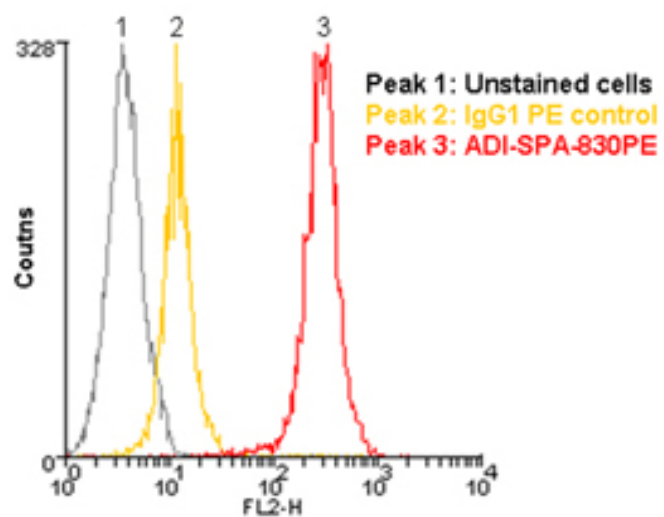
Ordering Information

[Order Online »](#)

ADI-SPA-830PE-D	50µg
ADI-SPA-830PE-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of 10^6 Jurkat cells stained using HSP90 mAb (AC88), R-Phycoerythrin Conjugate at a concentration of $10\mu\text{g/ml}$

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	Heat shock protein 90
Application	Flow Cytometry
Clone	AC88
Formulation	Liquid. In PBS containing 0.09% sodium azide.
Host	Mouse
Immunogen	Achlya ambisexualis (water mold) Hsp90.
Isotype	IgG1
Purity Detail	Protein G 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 ascites
Species Reactivity	Beluga, C. elegans, Chicken, Dog, E. coli, Fish, Fusobacterium nucleatum, Guinea pig, Hamster, Human, Monkey, Mouse, Mussel, Mycobacterium leprae, Porcine, Rabbit, Rat, Scallop, Sheep, Water mold
UniProt ID	Q8LLI6

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)