

HSP90 monoclonal antibody (AC88) (DyLight™ 488 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.

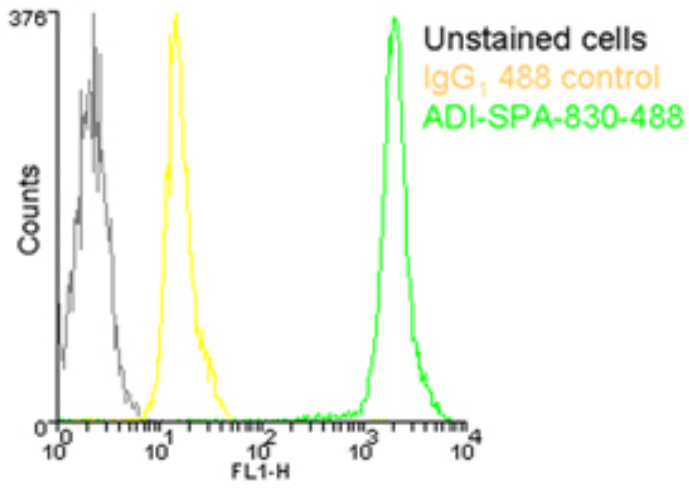
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 4

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of 10^6 Jurkat cells stained using HSP90 mAb (AC88), DyLight™ 488 Conjugate at a concentration of $50\mu\text{g/ml}$

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 (1:10) 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](#)

Last modified: May 29, 2024



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)