

HSP90 monoclonal antibody (2D11B9)

The 90kDa molecular chaperone family includes 90kDa heat shock protein Hsp90 and 94kDa glucose-regulated protein grp94, both major molecular chaperones of the cytosol and the endoplasmic reticulum. Mammalian cells contain isoforms Hsp90 α and Hsp90 β , encoded by separate genes. The amino acid sequences of human and yeast Hsp90 α are 85% and 90% homologous to those of Hsp90 β , 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 proteins 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 signalling and cell cycle control. In this context, researchers identified key regulatory proteins as substrates of Hsp90, including steroid receptors, cell cycle kinases involved in signal transduction, and p53. Hsp90 may act as a capacitor for morphological evolution by buffering widespread variation, potentially affecting morphogenic pathways. When temperature and other stress factors compromise *Drosophila* Hsp90 buffering, cryptic variant expression occurs, and selection can lead to the continued expression of these traits, even after Hsp90 function is restored.

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

Citations: 14

[View Online »](#)

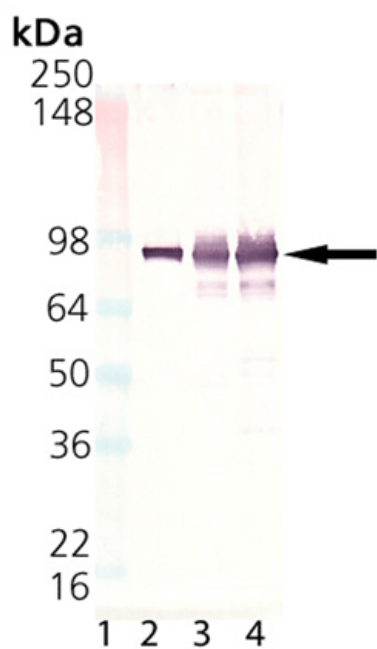
Ordering Information

[Order Online »](#)

ADI-SPA-831-050	50 μ g
ADI-SPA-831-200	200 μ g

Manuals, SDS & CofA

[View Online »](#)



Western Blot Analysis of HSP90, mAb (2D11B9) (Prod. No. ADI-SPA-831). Lane 1: MWM; Lane 2: HSP90 (human), (native) (Prod. No. ADI-SPP-770); Lane 3: HeLa, (cell lysate) (Prod. No. ADI-LYC-HL100); and Lane 4: HeLa (heat shocked), (cell lysate) (Prod. No. ADI-LYC-HL101).

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Heat shock protein 90

WB

Detects a band of ~92kDa by Western blot.

2D11B9

Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

Mouse

Recombinant human Hsp90.

IgG1

Protein G affinity purified.

Western Blot (1µg/ml, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Purified from ascites.

Dog, Drosophila, Fish, Guinea pig, Hamster, Horse,
Human, Monkey, Mouse, Porcine, Rat, Sheep

P07900 (HSP90 α), P08238 (HSP90 β)

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