

# HSP90α monoclonal antibody (9D2) (HRP 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 contain isoforms Hsp90 alpha and Hsp90 beta, 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 proteins in most cells. Hsp90 functions as part of the cell

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

Citations: 1

[View Online »](#)

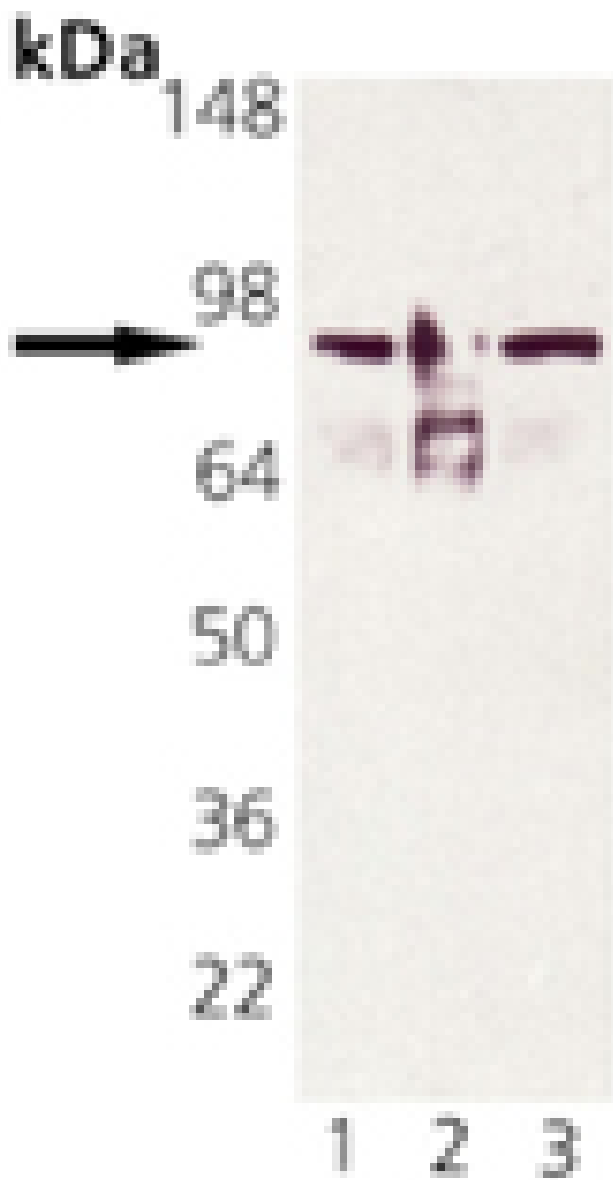
## Ordering Information

[Order Online »](#)

ADI-SPA-840HRP-D	50µg
ADI-SPA-840HRP-F	200µg

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of HSP90alpha mAb-HRP: Lane 1: HSP90 Native Protein, Lane 2: HSP90alpha Recombinant Protein, Lane 3: HeLa Cell Lysate (Heat Shocked).

## 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** HSP86, Heat shock protein 90α

**Application** WB

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

**Clone** 9D2

**Formulation** Liquid. In PBS, pH 7.2, containing 0.09% sodium azide diluted in proprietary stabilizer solution.

**Host** Rat

**Immunogen** Purified Hsp90 isolated from human therapeutic orchiectomy specimens.

**Isotype** IgG2a

**Purity Detail** Protein G affinity purified.

**Recommendation** Western Blot (1:2,000-1:50,000) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

**Source** Purified from rat ascites.

**Species Reactivity** Chicken, Human

**UniProt ID** P07900



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

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

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

France  
Phone: +33 1 42 44 55 55  
[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)

**Worry Free Guarantee** This antibody is covered by our **Worry Free Guarantee**