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 »

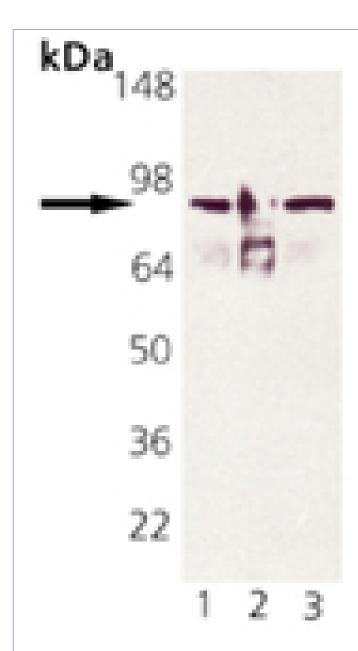
Ordering Information

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 NotesDetects 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.

lgG2a

Purity Detail Protein G affinity purified.

Recommendation Dilutions/ConditionsWestern 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

_

Last modified: August 6, 2025

