

HSP90β (human), (recombinant)

Recombinant human Hsp90β, a constitutively expressed cytosolic chaperone that stabilizes and folds key signaling proteins, steroid receptors, and

kinases, supporting cell cycle control and stress responses in an ATP-dependent manner.

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 express inducible HSP90α and constitutive HSP90β isoforms that are encoded by separate genes. The amino acid sequences of human and yeast HSP90α are 85% and 90% homologous to those of HSP90β, 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 signaling and cell cycle control by serving as a chaperone for many key signaling molecules including steroid receptors, cell cycle kinases involved in signal transduction, and p53. As many of these client proteins are known oncogenes, Hsp90 inhibitors such as 17-AAG, a geldanamycin analog, have been of benefit in the treatment of many cancers.

Citations: 7

[View Online »](#)

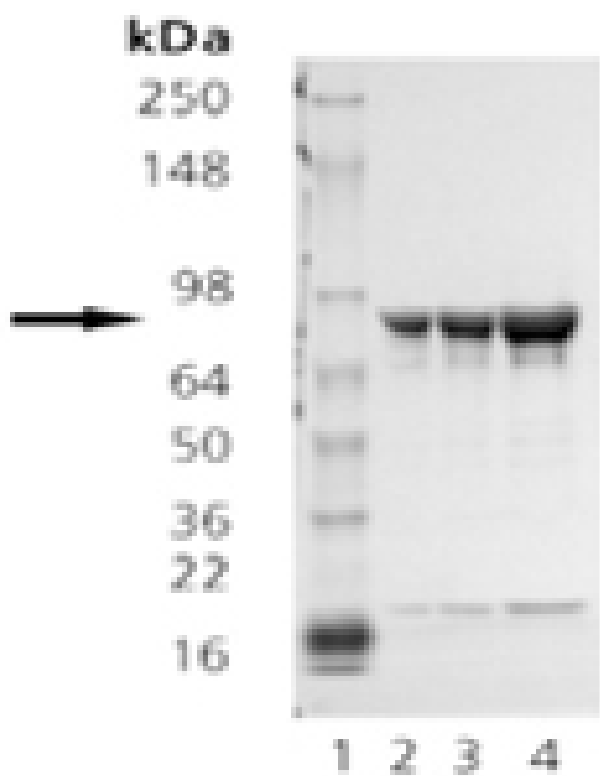
Ordering Information

[Order Online »](#)

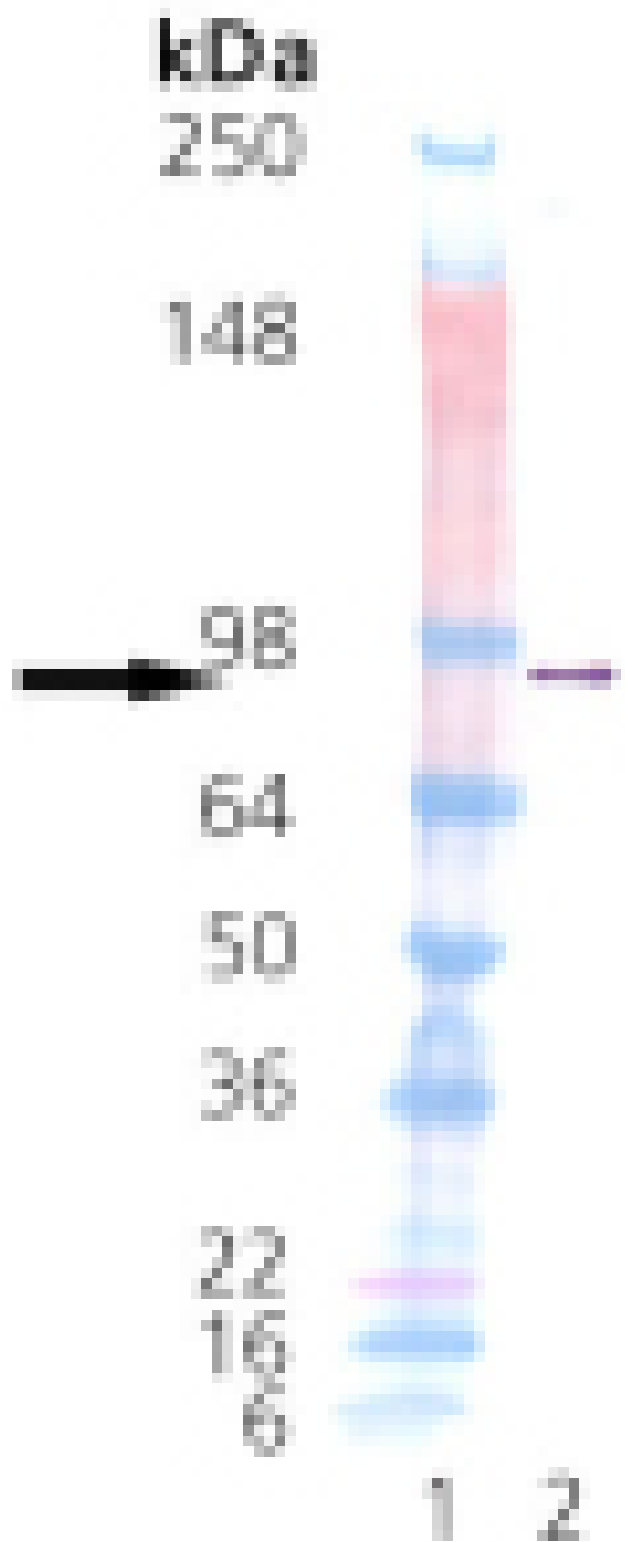
ADI-SPP-777-D	50μg
ADI-SPP-777-F	200μg

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis: Lane 1: MW marker, Lane 2: 1 μ g, Lane 3: 2 μ g, Lane 4: 5 μ g HSP90b



Western Blot analysis: Lane 1: MWM, Lane 2: HSP90 beta recombinant protein probed with HSP90 beta, mAb.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name HSP84, Heat shock protein 90β

Application Notes Western blot control.

Formulation Liquid. In 20mM TRIS, pH 7.5, containing 500mM sodium chloride, 5% glycerol, and 1mM DTT.

MW ~90kDa

Purity ≥85% (SDS-PAGE; Western blot)

Purity Detail Purified by multi-step chromatography.

Source Produced in *E. coli*.

UniProt ID P08238

Last modified: October 9, 2025



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)