HSP90 (human), (native)

Native human Hsp90, a highly conserved cytosolic chaperone that, in cooperation with co-chaperones, stabilizes client proteins including kinases, transcription factors, and

The Hsp90 camely express shock proteins segressed as one of the most abundantly expressed and highly conserved families of cellular chaperones dependent folding and activation. 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, bcrabl, p53, and hTERT, are members of well characterized oncogenic pathways, making Hsp90 inhibitors useful anticancer agents.

Citations: 13

View Online »

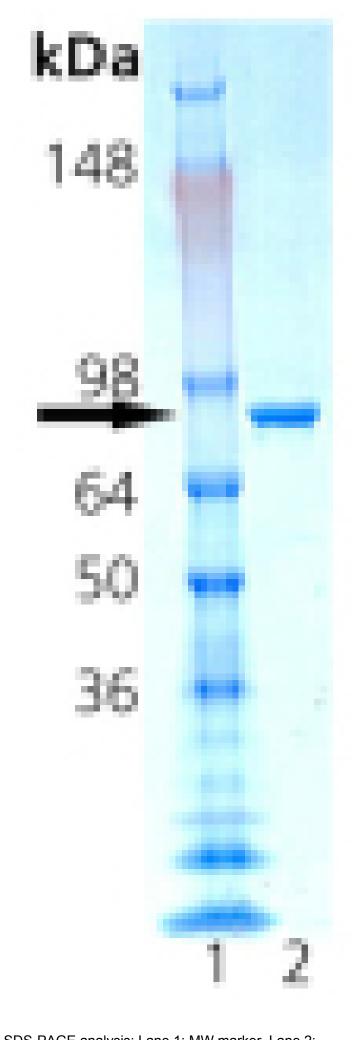
Ordering Information

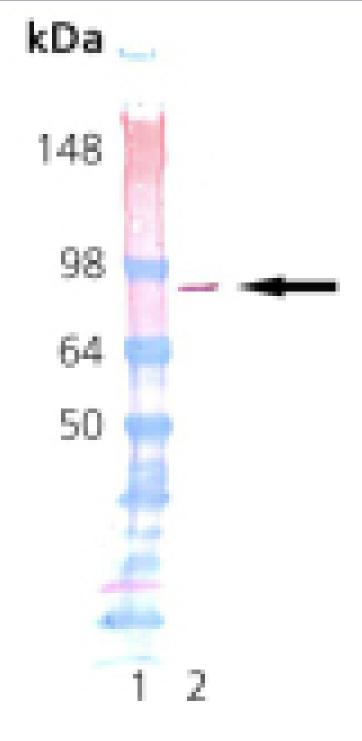
Order Online »

ADI-SPP-770-D	50µg
ADI-SPP-770-F	200µg

Manuals, SDS & CofA

View Online »





Western Blot analysis: Lane 1: MWM, Lane 2: Natural Source Human HSP90 Protein.

SDS-PAGE analysis: Lane 1: MW marker, Lane 2: 1.0µg of Natural Source Human HSP90 Protein.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name HSP86, Heat shock protein 90

Application Notes Western blot control.

Formulation Liquid. In 50mM TRIS-HCl, pH 7.5, containing 1.0mM DTT

and 100mM sodium chloride.

MW ~90kDa

Purity ≥90% (SDS-PAGE; Western blot)

Purity Detail Purified by multi-step chromatography.

Source Produced in HeLa cells.

UniProt ID P07900 (HSP90α), P08238 (HSP90β)

Last modified: October 9, 2025



eu@enzolifesciences.com