

# HSP90 $\alpha$ (human), ELISA kit

**Most sensitive, rapid ELISA kit for highly specific, quantitative detection of human HSP90 $\alpha$ .**

The HSP90 $\alpha$  (human), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in < 3 hours. Absorbance is read at 450 nm. Save time, money, and precious sample with fully quantitative results and increased sensitivity compared to Western blot analysis.

The Hsp90 family of heat shock proteins represents one of the most abundantly expressed and highly conserved families of cellular chaperones 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, bcr-abl, p53, and hTERT, are members of well characterized oncogenic pathways, making Hsp90 inhibitors useful anticancer agents.

Citations: 29

[View Online »](#)

## Ordering Information

[Order Online »](#)

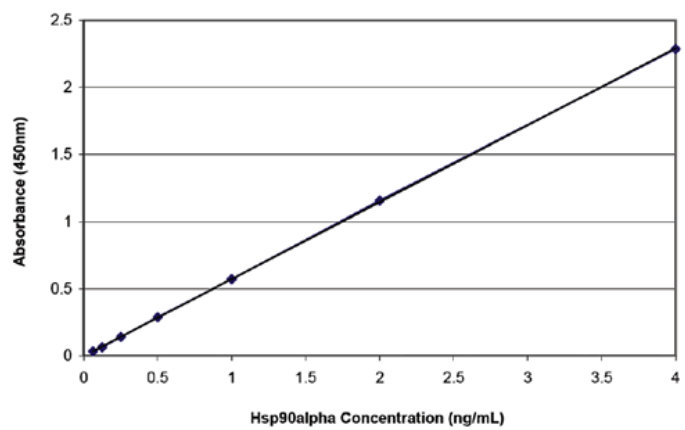
ADI-EKS-895

96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Reproducible with less than 10% variation between assays
- Highly sensitive detecting as little as 50 pg/ml of human HSP90 $\alpha$
- Higher throughput format, test up to 40 samples in duplicate in just 3 hours
- Fully quantitative results that surpass semi-quantitative Western blot analysis



## Handling & Storage

**Use/Stability** Store all components at 4°C, except standard at -20°C.

**Shipping** Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Heat shock protein 90α, HSP86

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of HSP90α in cell lysates, serum, and tissue of human origin.

**Assay Time** <3 hours

**Compatibility** This product is compatible with the [Absorbance 96 Plate Reader](#).

**Contents** Microtiter plate, 5X Extraction reagent 2, Standard, Sample diluent, Wash buffer concentrate, HRP Conjugate, Conjugate diluent, TMB Substrate, Stop solution 2

**Crossreactivity** HSP90α (100%). No reactivity with HSP90β, Grp94, HSP60, or HSP70 (HSP72).

**Sensitivity** 50 pg/ml (range 62.5 – 4,000 pg/ml)

**Species Reactivity** Human

**UniProt ID** P07900

**Wavelength** 450 nm

**Last modified: May 29, 2024**



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-be@enzolifesciences.com)

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