

HSP27 (human), ELISA kit

Most comprehensive portfolio of heat shock protein (chaperone) products commercially available.

The HSP27 (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. Highly sensitive quantitative determination of Hsp27 independent of phosphorylation state

Hsp27 is one of the most common members of the highly conserved and ubiquitously expressed family of small heat shock proteins (sHsp), which also includes alphaB-crystallin. It is characterized by a conserved C-terminal alpha-crystallin domain consisting of two anti-parallel beta-sheets that promote oligomer formation required for its primary chaperone function as inhibitor of irreversible protein aggregation. Hsp27 oligomerization is modulated by post-translational phosphorylation of Hsp27 at three serine residues, Ser15, Ser78, and Ser82, by a variety of protein kinases including MAPKAPK-3, PKAc- α , p70 S6K, PKD I, and PKC- δ . Hsp27 has been shown to inhibit actin polymerization by binding of unphosphorylated Hsp27 monomers to actin intermediate filaments. Anti-apoptotic functions of Hsp27 have also been identified through interactions with DAXX7, activation of Akt, and inhibition of apoptosome formation. Evidence suggests altered expression of Hsp27 is implicated in the pathogenesis of breast, ovarian, and prostate cancer.

Citations: 21

[View Online »](#)

Ordering Information

[Order Online »](#)

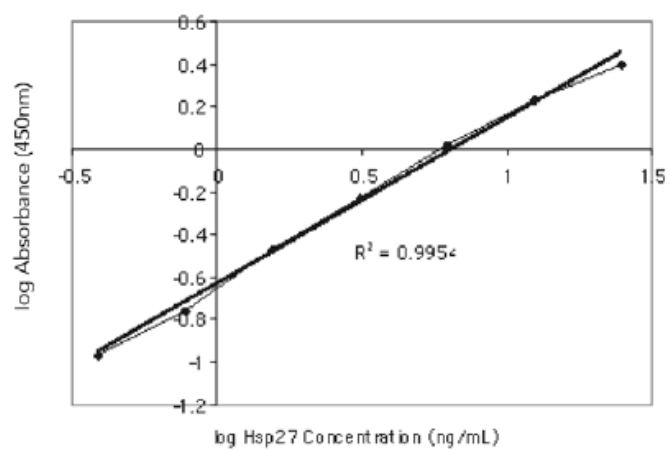
ADI-EKS-500

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Highly sensitive ELISA, measuring as little as 0.39 ng/ml
- Fast results from up to 39 samples in duplicate in less than 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 27

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of human HSP27 in cell lysates, plasma, serum, and tissue.

Assay Time <3 hours

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

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

Crossreactivity HSP27 (100%). No reactivity with HSP25 (mouse or rat) or α B-Crystallin.

GenBank ID L39370

Sensitivity 0.39 ng/ml (range 0.39 – 25 ng/ml)

Species Reactivity Human

UniProt ID P04792

Wavelength 450 nm



ENZO LIFE SCIENCES,
INC.
Phone: +1 949 946 9120
info@enzolifesciences.com
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info@enzolifesciences.com
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com
be@enzolifesciences.com

France
Phone: +33 472 440 655
info@enzolifesciences.com
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com
de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info@enzolifesciences.com
uk@enzolifesciences.com