

IMMUNOSET[®] HSP27

high sensitivity (human), ELISA development set

The IMMUNOSET[®] HSP27 high sensitivity (human) EIA development set contains sufficient reagents for the development of 5 x 96 well EIA plates which provides analysis of ~200 samples in duplicate. Save time, money, and sample by developing your own EIA kits. Simply coat your plates and go! This kit provides superior accuracy with fully quantitative results compared to Western blot analysis and has low reactivity with related molecules.

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 α B-crystallin. It is characterized by a conserved C-terminal α -crystallin domain consisting of two anti-parallel β -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: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

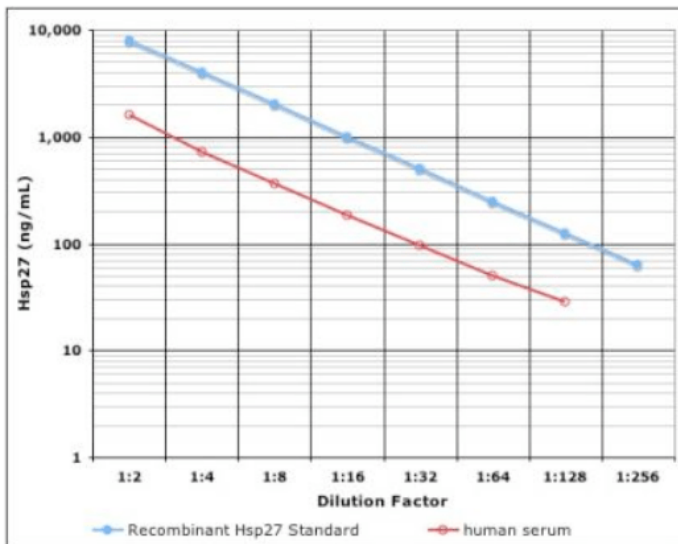
ADI-960-076

5x96 wells

Manuals, SDS & CofA

[View Online »](#)

- Cost effective - coat your own plates and stretch your research budget
- Quantitative - obtain fully quantitative statistically significant results
- Ultra-Sensitive - detect as little as 97 pg/mL Hsp27 compared to microgram levels in Western blot
- Reliable - obtain accurate results with thoroughly tested reagents in multiple complex sample matrices



Handling & Storage

Long Term Storage +4°C

Shipping Blue 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, and serum.
Contents	Capture antibody, Standard, Detection antibody, SA-HRP
Crossreactivity	Specific for HSP27. Does not cross-react with HSP25, Cpn10, HO-1, HSP40, or αB-crystallin.
GenBank ID	L39370
Sensitivity	97 pg/ml (range 0.1 – 3.2 ng/ml)
UniProt ID	P04792
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-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)