

# [pSer<sup>15</sup>]HSP27 (human), ELISA kit

**Only commercially available ELISA kit for ultra-sensitive detection of human phospho-Ser15 HSP27 in cell lysates, plasma, and serum.**

The [pSer<sup>15</sup>]HSP27 (human), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in 3 hours. Save time, money, and precious sample with fully quantitative results and increased sensitivity compared to Western blot analysis.

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- $\alpha$ , p70 S6K, PKD I, and PKC- $\delta$ . 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.

## Ordering Information

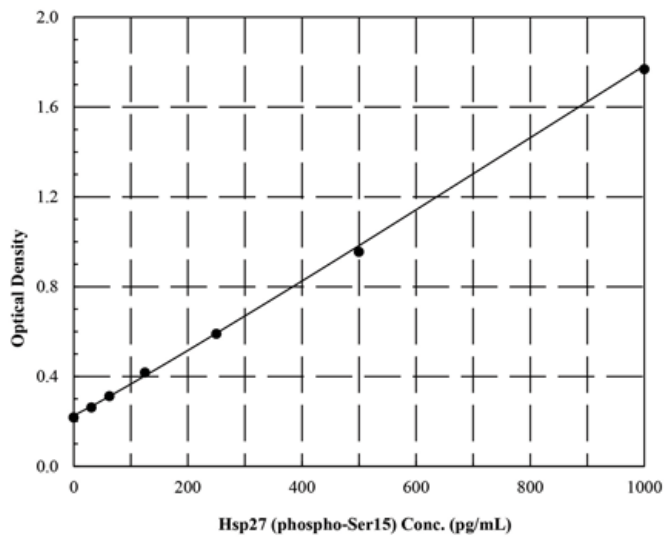
[Order Online »](#)

ADI-900-170	96 wells
-------------	----------

## Manuals, SDS & CofA

[View Online »](#)

- Ultra-sensitive and specific - be confident with your results and use less sample per test
- Quantitative - obtain statistically significant results compared to semi-quantitative Western blot analysis
- Higher throughput - assay up to 40 samples in duplicate in just 3 hours



# Handling & Storage

**Use/Stability** Store all components at 4°C, except standard and antibody at -20°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 [pSer<sup>15</sup>]  
]HSP27 in cell lysates, plasma, and serum.

**Assay Time** 3 hours

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

**Contents** Microtiter plate, Conjugate, Antibody, Assay buffer 27, Wash buffer concentrate, Standard, TMB Substrate, 5X Extraction reagent, Stop solution 2

**Crossreactivity** Phospho HSP27 (Ser15) (100%), HSP27 (non-phospho) (0.885%), HSP70 (0.023%) and <0.016%: HSP25, HSP40, HSP90, αB Crystallin, αA Crystallin

**GenBank ID** L39370

**Sensitivity** 10.15 pg/ml (range 31.25 – 1000 pg/ml)

**Species Reactivity** Human

**UniProt ID** P04792

**Wavelength** 450 nm



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.948.0120  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

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