

[pSer78]HSP27 (human), ELISA kit

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

The [pSer⁷⁸]HSP27 (human), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in 3 hours.

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 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: 3 [View Online »](#)

Ordering Information [Order Online »](#)

ADI-900-165	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 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 [pSer⁷⁸]
]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 (Ser78) (100%), HSP27 (non-phospho) (0.1%) and <0.03%: HSP25, HSP40, HSP70, HSP90, αB Crystallin, αA Crystallin

GenBank ID L39370

Sensitivity 4.30 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
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