

HSP27 (phospho) (human), (recombinant)

Recombinant human phospho-Hsp27 that functions as a small heat shock protein, with

phosphorylation-modulated oligomerization, preventing protein aggregation, regulating cytoskeletal dynamics, and providing anti-apoptotic protection.

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 antiparallel β -sheets that promote oligomer formation required for its primary chaperone function as inhibitors of irreversible protein aggregation. Hsp27 oligomerization is modulated by post-translational phosphorylation of Hsp27 at three serine residues, Ser-15, Ser-78, and Ser-82, by a variety of protein kinases including MAPKAPK-3, PKA α , p70S6K, PKD I, and PKC δ . Hsp27 has been shown to inhibit actin polymerization by binding of unphosphorylated Hsp27 monomers to actin intermediate filaments. Antiapoptotic 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: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-SPP-716-F	200 μ g
---------------	-------------

Manuals, SDS & CofA

[View Online »](#)

kDa

148

98

64

50

36

22

16

6

1

2



kDa

250

148

98

64

50

36



22

16

6

1

2

3

4

5

SDS-PAGE analysis: Lane 1: MW marker, Lane 2: 0.5µg, Lane 3: 1µg, Lane 4: 2µg, Lane 5: 5µg Hsp27 (phospho) protein

Western Blot analysis: Lane 1: MWM, Lane 2: Hsp27 (phospho) protein, probed with Hsp27 (pSer82) pAb

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name HspB1, Heat shock protein 27

Application Notes Western blot control.

Formulation Liquid. In 60mM Tris-HCl, pH 7.5, containing 5mM sodium chloride, 1mM EGTA, 0.5mM EDTA, 15mM DTT, 0.1mM ATP, and 15mM magnesium chloride.

GenBank ID L39370

MW ~27kDa

Purity ≥90% (SDS-PAGE; Western blot)

Purity Detail Purified by multi-step chromatography.

Source Produced in *E. coli*.

UniProt ID P04792

Last modified: October 9, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

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

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

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