

# [pSer<sup>15</sup>]HSP27 polyclonal antibody

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.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 4

[View Online »](#)

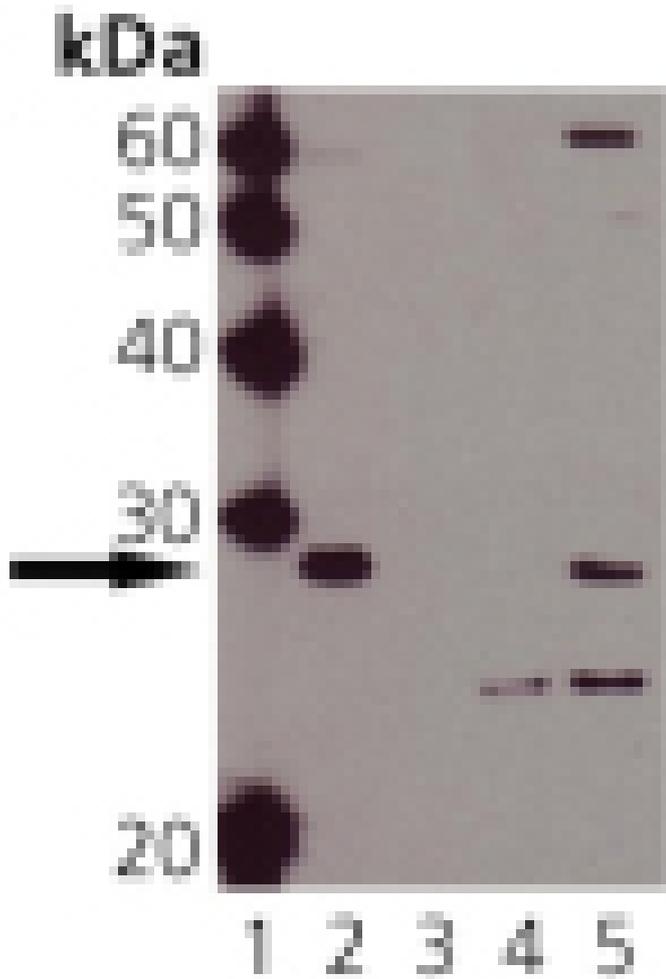
Ordering Information

[Order Online »](#)

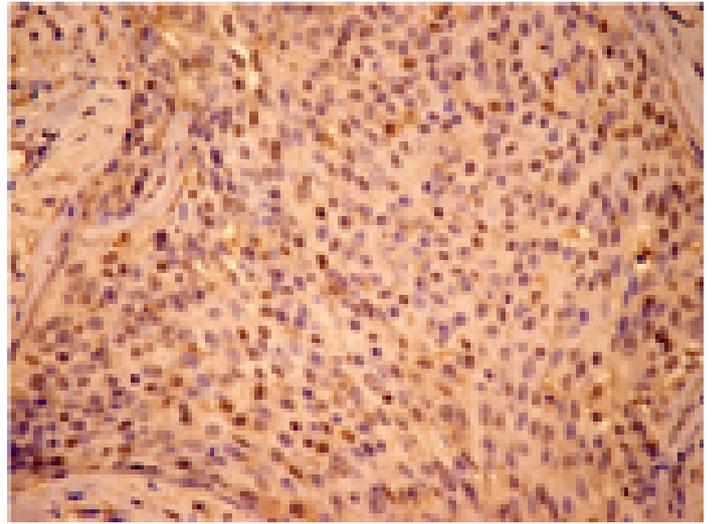
ADI-SPA-525-D	50µg
---------------	------

Manuals, SDS & CofA

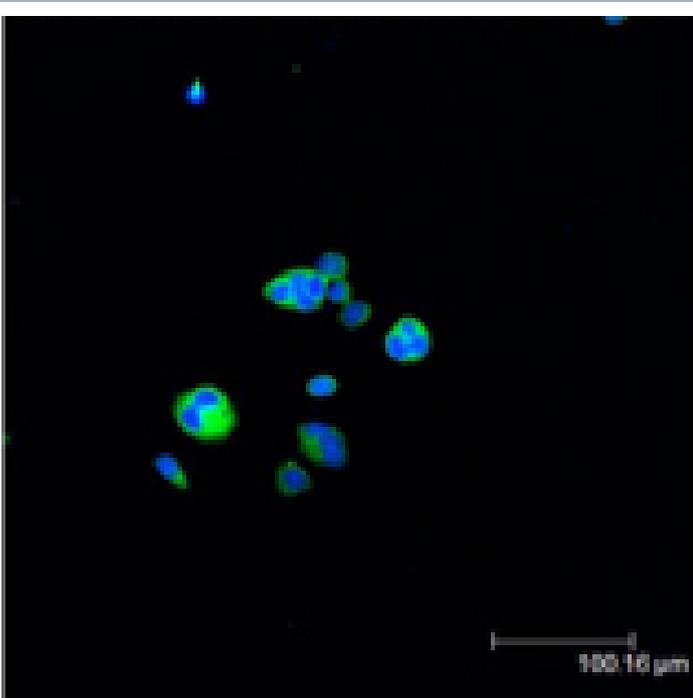
[View Online »](#)



Western blot analysis of HSP27 (pSer15): Lane 1: MW marker, Lane 2: HSP27 (phospho) Recombinant Human Protein, Lane 3: HSP27 Recombinant Human Protein, Lane 4: HeLa, Lane 5: HeLa, Heat-Shocked.



Immunohistochemistry analysis of human breast cancer tissue with Hsp27 (pSer15) pAb



Immunofluorescent analysis (confocal) staining of HeLa cells using Hsp27 (pSer15) pAb (green); nuclei are stained in blue pseudocolor using DRAQ.

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** HspB1, Heat shock protein 27

**Application** IF, IHC, WB

**Application Notes** Detects a band of ~27kDa by Western blot.

**Formulation** Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

**GenBank ID** L39370

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to a portion of human HSP27 phosphorylated at Ser15.

**Purity Detail** Peptide affinity purified.

**Recommendation Dilutions/Conditions** Immunohistochemistry (1:50)Western Blot (1:1,000, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Source** Purified from rabbit serum.

**Species Reactivity** Bovine, Dog, Guinea pig, Human, Monkey

**UniProt ID** P04792

**Worry-free Guarantee** This antibody is covered by our [Worry-Free Guarantee](#)

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)