

[pSer⁸²]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: 5

[View Online »](#)

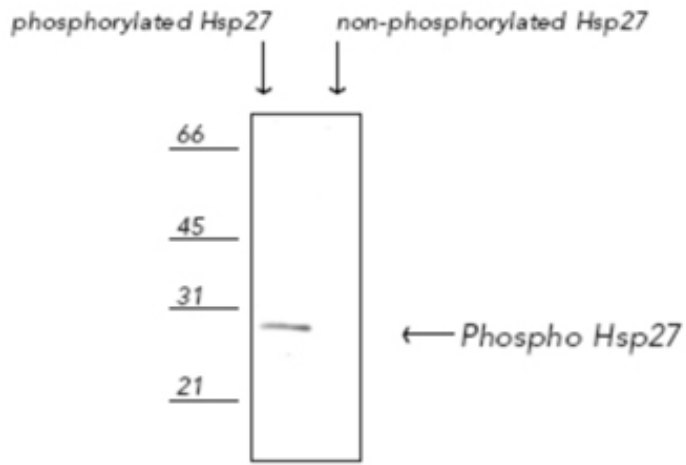
Ordering Information

[Order Online »](#)

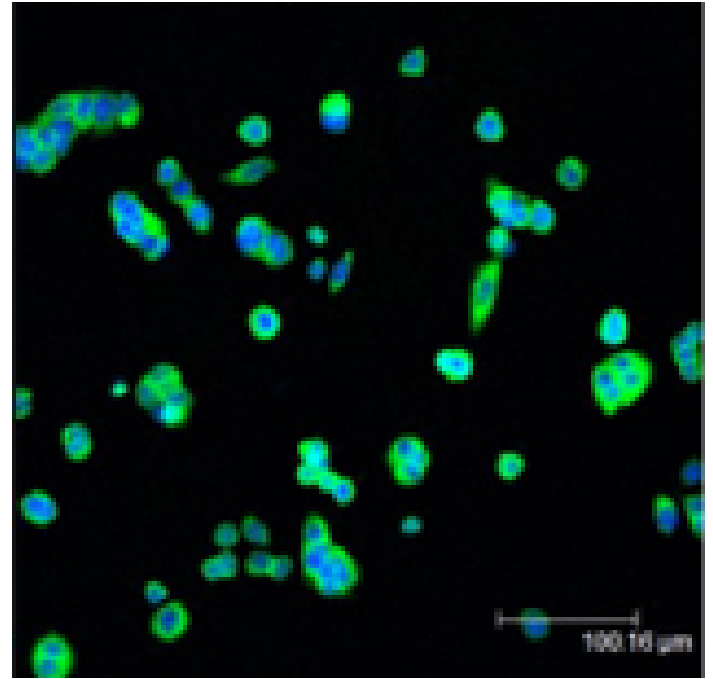
ADI-SPA-524-D	50µl
---------------	------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of phosphorylated HSP27 and nonphosphorylated HSP27 proteins probed with HSP27 (pSer82) pAb



Immunofluorescent analysis (confocal) staining of HeLa cells using Hsp27 (pSer82) pAb (green); nuclei are stained in blue pseudocolor using DRAQ5

Handling & Storage

Use/Stability	Serum should be stored at -20°C for long term storage. For short term storage, 4°C, supplement with 0.1% sodium azide.
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, IP, WB
Application Notes	Detects a band of ~27kDa by Western blot.
Formulation	Liquid. Whole rabbit serum.
GenBank ID	L39370
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of human HSP27 phosphorylated at Ser82.
Recommendation Dilutions/Conditions	Immunoprecipitation (1:100)Western Blot (1:2,000, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Whole rabbit serum.
Species Reactivity	Bovine, Chicken, Dog, Guinea pig, Human, Mouse, Rat, Sheep
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)