[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 »

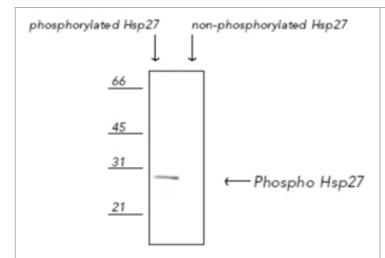
Ordering Information

Order Online »

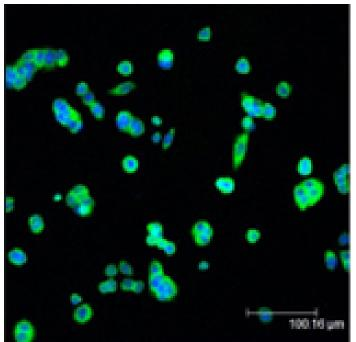
ADI-SPA-524-D	50μΙ
ADI-SPA-524-F	200μΙ

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 -20oC for long term storage. For short term storage, 4oC,

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 NotesDetects 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 ReactivityBovine, 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



info-

eu@enzolifesciences.com