HSP27 monoclonal antibody (G3.1) (biotin conjugate)

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: 1

View Online »

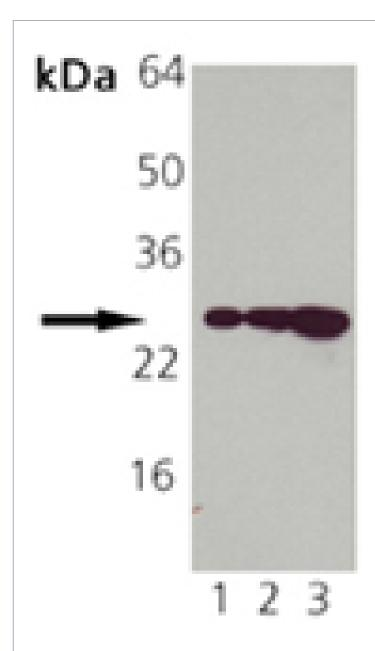
Ordering Information

Order Online »

ADI-SPA-800B-D	50µg
ADI-SPA-800B-F	200μg

Manuals, SDS & CofA

View Online »



Western blot analysis of Biotinylated HSP27 mAb: Lane 1: HSP27 Recombinant Protein, Lane 2: HeLa Cell Lysate (Heat Shocked), Lane 3: Vero (Heat Shocked).

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 WB

Application NotesDetects a band of ~27kDa by Western blot.

Clone G3.1

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09%

sodium azide.

GenBank ID L39370

Host Mouse

Immunogen Native human HSP27.

lsotype IgG1

Purity Detail Protein G affinity purified.

Recommendation Dilutions/ConditionsWestern 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 mouse ascites.

Species Reactivity Bovine, Fish, Human, Monkey, Mouse, Rat

UniProt ID P04792

Last modified: May 29, 2024



info-

eu@enzolifesciences.com