HSP27 (phospho) (human), (recombinant)

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

pho25 is non-of the most-common members of the highly conserved and ubiquitously expressed family of small heat shock proteins (sHSP), which proventing protein aggregation, regulating also includes aB-crystallin. It is characterized by a conserved C-terminal acrystalline consisting io swall parallel is need that promote eligenetic matter promote eligenetic matter 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, PKAcα, p70S6K, PKD I, and PKCδ. 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 »

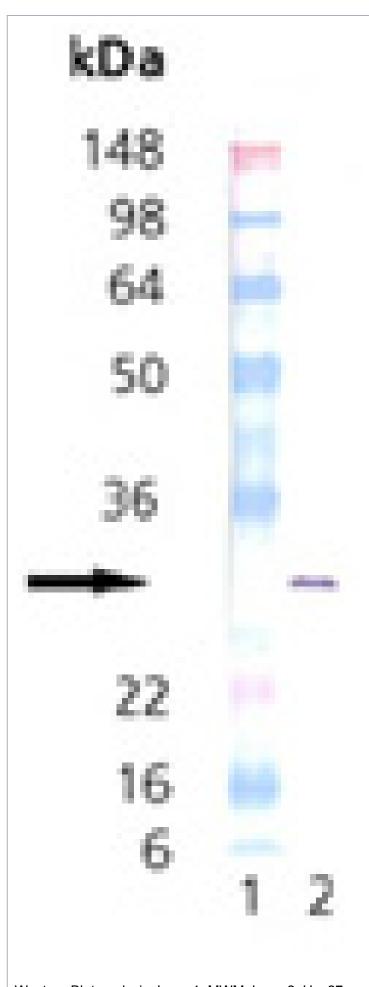
Ordering Information

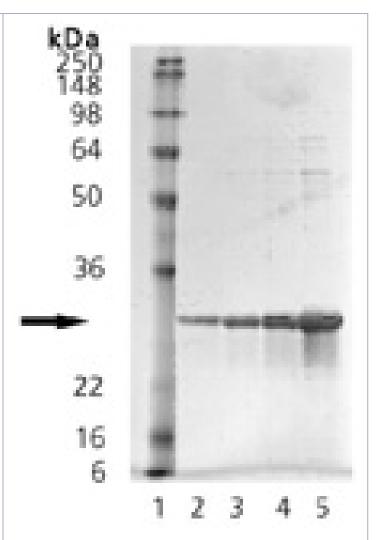
Order Online »

ADI-SPP-716-D	50µg
ADI-SPP-716-F	200μg

Manuals, SDS & CofA

View Online »





SDS-PAGE analysis: Lane 1: MW marker, Lane 2: 0.5 μ g, Lane 3: 1 μ g, Lane 4: 2 μ g, Lane 5: 5 μ g 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

