HSP27 (human), (recombinant)

Recombinant human Hsp27 that functions as a small heat shock protein, preventing protein aggregation, supporting refolding with Hsp70, regulating cytoskeletal

thynan Hsp27, another Hsp25 and ab constallin are past of a diverse family of small heat shock proteins which are produced in all organisms. They protection. In the protein aggregation of the protein aggregation of the protein aggregation. Hsp27 is believed to exist mainly as oligomers of as many as 8-40 Hsp27 protein monomers in cells. Data suggests that the large oligomers of Hsp27 have a chaperone-like activity by serving as a site where unfolded proteins may bind until ATP and Hsp70-dependent refolding can occur. The phosphorylation and oligomerization state of Hsp27 has been suggested to regulate microfilament organization, since only the nonphosphorylated lower molecular weight forms of Hsp27 bind actin barbed ends and inhibit polymerization. Hsp27 is also believed to protect cells by enhancing cellular glutathione levels, as elevated glutathione levels have been measured in cells overexpressing Hsp27. Data from studies using wild-type Hsp27 and mutant forms in which the serine phosphorylation sites were

Citations: 15

View Online »

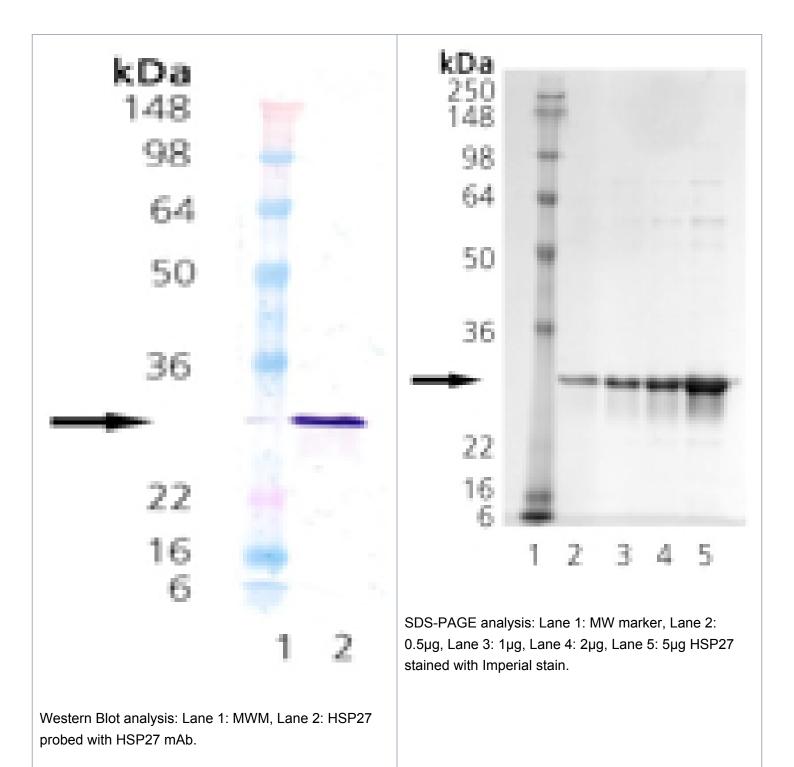
Ordering Information

Order Online »

ADI-SPP-715-D	50µg
ADI-SPP-715-F	200μg

Manuals, SDS & CofA

View Online »



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 20mM TRIS, pH 7.5, containing 10mM sodium

chloride, 1mM EDTA, and 1mM DTT.

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



eu@enzolifesciences.com