

# HSP27 (human), (recombinant)

Human Hsp27, mouse Hsp25 and ab-crystallin are part of a diverse family of small heat shock proteins which are produced in all organisms. They function as chaperone-like proteins by binding unfolded polypeptides and preventing uncontrolled 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 »](#)

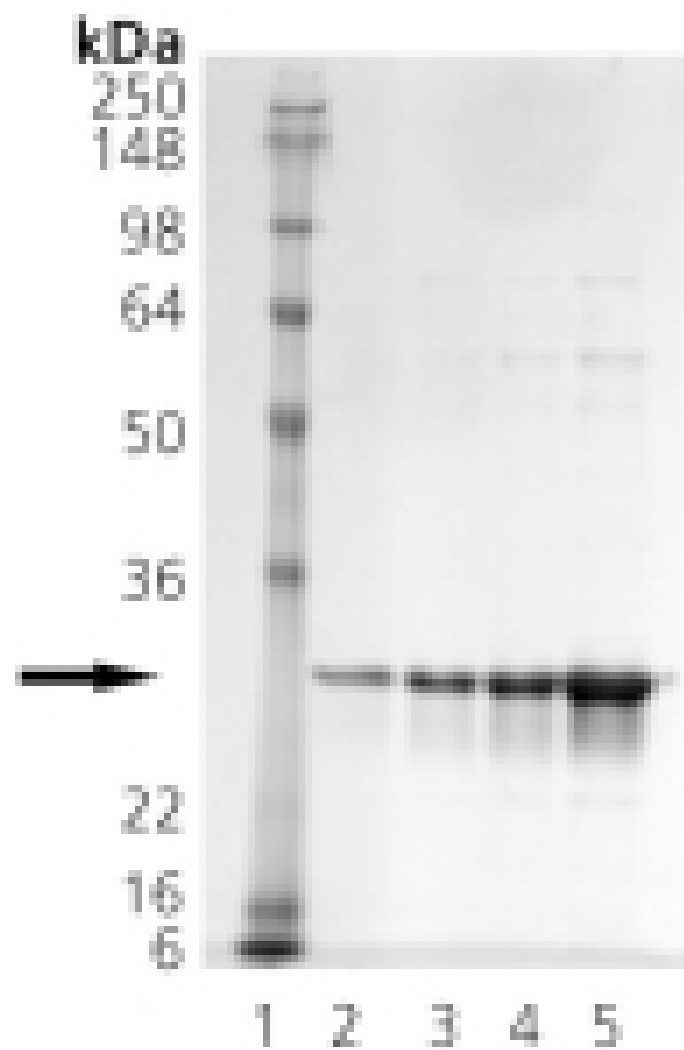
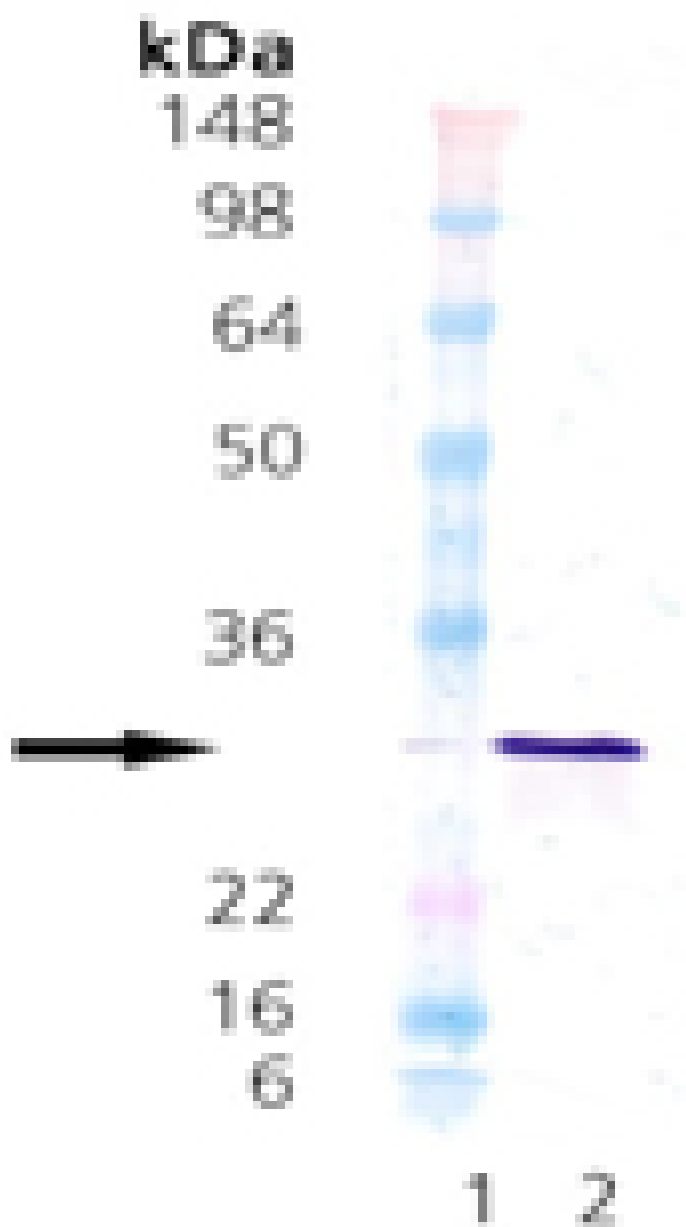
## Ordering Information

[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



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)