

HSP27 (low endotoxin) (human), (recombinant)

Recombinant human HSP27 (low endotoxin) that functions as a molecular chaperone and inhibitor of protein aggregation.

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.

Citations: 2

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

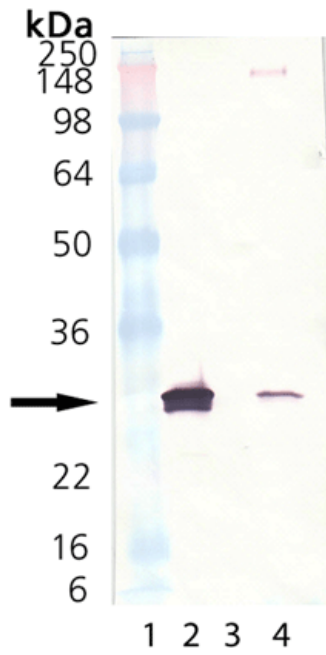


Figure 2: Western Blot Analysis – Lane 1: MW Marker; Lane 2: 100ng Hsp27; Lane 3: 20 μ g HeLa lysate; Lane 4: 20 μ g Heat-shocked HeLa lysate.

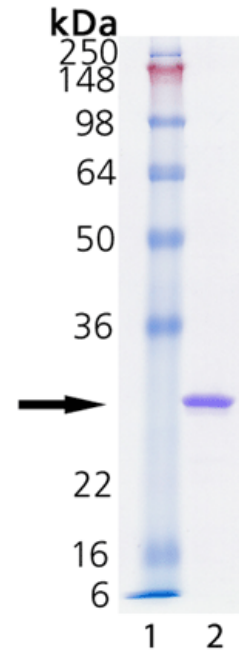


Figure 1: SDS-PAGE Analysis – Lane 1: MW Marker; Lane 2: 1 μ g Hsp27

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. |
| Endotoxin Content | <50EU/mg purified protein (LAL test) |
| Formulation | Liquid. In Dulbecco's PBS. |
| GenBank ID | L39370 |
| MW | ~27kDa |
| Purity | ≥90% (SDS-PAGE; Western blot) |
| Purity Detail | Purified by multi-step chromatography. |
| Source | Produced in <i>E. coli</i> . |
| UniProt ID | P04792 |

Last modified: October 9, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com