

HSP25 (mouse), (recombinant)

Recombinant mouse HSP25 that acts as a small heat shock protein, preventing protein aggregation and supporting cellular stress responses.

Mouse Hsp25, human Hsp27, and $\alpha\beta$ -crystallin are part of a diverse family of small heat shock proteins (sHsps) which are produced in all organisms. The overall homology between the different sHsps is low and they are grouped together based on conserved sequences in the C-terminal half of the protein and short conserved phenylalanine-rich stretches near the N terminus. Mammalian sHsps are expressed constitutively under physiological conditions but stress factors such as heat shock induce a strong up-regulation of protein levels by 10-20-fold to maximum concentrations of 0.1% of the cellular protein. They function as chaperone-like proteins by binding unfolded polypeptides and preventing uncontrolled protein aggregation. sHsps all share the striking feature of forming high molecular weight oligomeric complexes of variable size. The quaternary structure of sHsps is essential for their function and regulation of activity but its basic properties are still rather poorly understood. Data indicates that Hsp25 is a dynamic tetramer of tetramers with a unique ability to refold and reassemble into its active quaternary structure after denaturation. Current studies demonstrate that Hsp25 helps facilitate the glutathione-redox cycle by enhancing glutathione utilization and maintaining the cellular glutathione pool in favor of the reduced states.

Citations: 3

[View Online »](#)

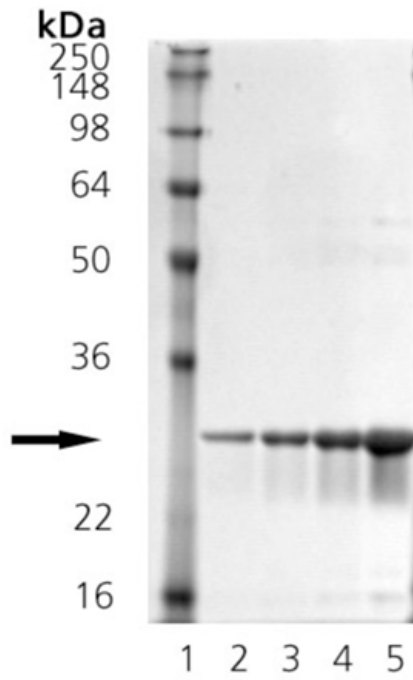
Ordering Information

[Order Online »](#)

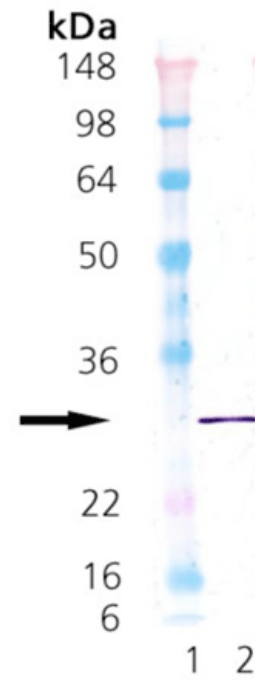
ADI-SPP-510-D	50 μ g
---------------	------------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis: Lane 1: MWM, Lane 2: 0.5 μg, Lane 3: 1 μg, Lane 4: 2 μg, Lane 5: 5 μg Hsp25.



Western Blot analysis: Lane 1: MW marker, Lane 2: HSP25 protein probed with anti-HSP25 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 25
Application Notes	Western blot control.
Formulation	Liquid. In Dulbecco's PBS.
GenBank ID	L07577
MW	~28kDa
Purity	≥90% (SDS-PAGE; Western blot)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> .
UniProt ID	P14602

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