

HSP25 polyclonal antibody

Hsp25 is a member of the family of small heat shock proteins (sHsps), an evolutionarily conserved family of stress proteins which exhibit molecular weights between 14-45 kDa. The sHsps are related to the alpha-crystallins, and generally exist as large oligomeric complexes that serve to stabilize and assist in refolding of denatured proteins. Hsp25 is highly homologous to the human Hsp27, and its expression is seen in many species, including mouse, rat, bovine, canine, and hamster cells.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 80

[View Online »](#)

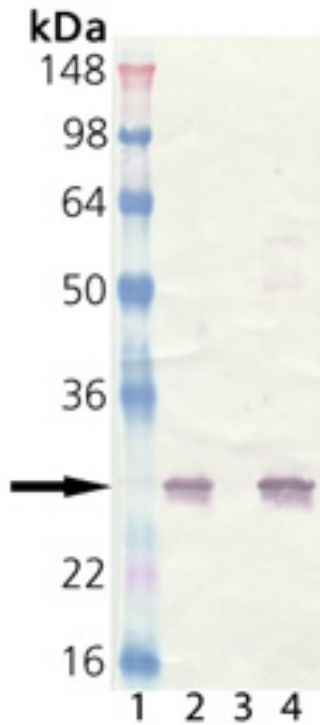
Ordering Information

[Order Online »](#)

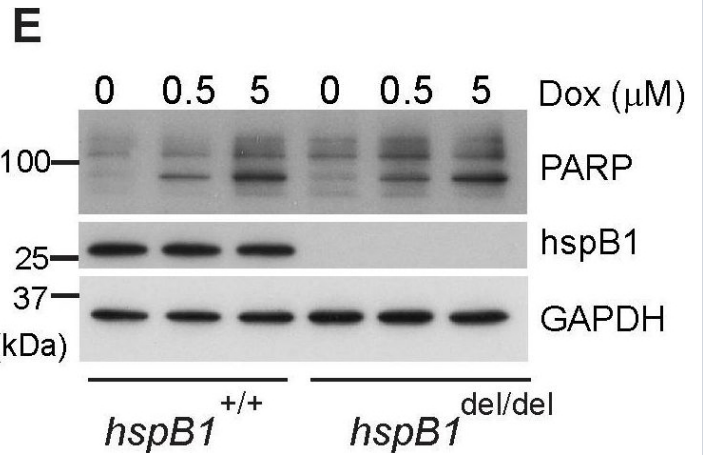
ADI-SPA-801-D	50µg
ADI-SPA-801-F	200µg

Manuals, SDS & CofA

[View Online »](#)



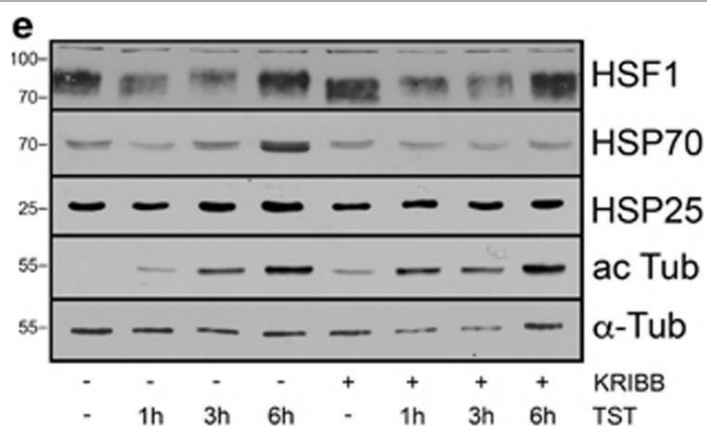
Western Blot analysis of HSP25, pAb. Lane 1: MW Marker; Lane 2: PC-12 Cell Lysate (Heat Shock), Lane 3: 3T3 Cell Lysate (Heat Shock), Lane 4: HSP25 (mouse), (recombinant) (Prod. No. ADI-SPP-510).



E

HspB1 deficiency increases IL-1-induced IL-6 expression and inhibits proliferation in fibroblasts. A, Primary wild-type and hspB1del/del MEF were treated with IL-1 (20 ng/ml) for 4 h or left untreated. Graph shows the concentration of IL-6 in culture medium as determined by ELISA and normalised against values for IL-1-treated wild-type MEF for three separate batches of cells (* $P < 0.05$). B, MEF were treated as in (A), lysed, RNA extracted and IL-6 and GAPDH mRNAs quantified by qRT-PCR. Plot shows IL-6 mRNA/GAPDH mRNA normalised to the value for wild-type cells treated with IL-1 for 1 h. C, Growth curve analysis (means \pm SEM) of wild-type and hspB1del/del MEF determined by MTT assay at different days post-seeding ($n = 3$; * $P < 0.05$, ** $P < 0.01$, *** $P < 0.001$) or by counting trypan-blue excluded cells ($n = 2$). D, Plot of mean (%) TUNEL-positive (\pm SEM) cells for three different batches of MEF per genotype. E, MEF were treated with different concentrations of doxorubicin (as indicated) for 8 h to induce apoptosis, or left untreated, cells lysed and lysates analysed by western blot for the full-length and the cleaved form of PARP. Similar results were obtained in three independent experiments.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Heat shock protein B1-deficient mice display impaired wound healing. *PLoS One* (2013)



Tubastatin A induces heat-shock protein expression by activating heat-shock factor 1. 661W cells were treated with 1, 5 and 10 μM of tubastatin A (TST) for 24 h (a), or in (b) with 10 μM TST for 8 h, or with 200 μM H₂O₂ for 6 h or were preincubated with 10 μM TST for 2 h followed by incubation with 200 μM H₂O₂ for 6 h. ac Tub, acetylated tubulin. α -Tub, α -tubulin. Co, untreated control. Quantitative evaluation of immunoblot analysis revealed a significant increase in heat-shock protein (HSP) 70 level after 8 h (c), while HSP25 was significantly enhanced after 24 h (d); n=4. (e) Heat-shock factor 1 (HSF1) activity was investigated using immunoblot analysis of 661W cell extracts that were treated 10 μM TST for 1, 3 and 6 h, or with 5 μM KRIBB11 (KR) for 6.5 h alone, or preincubated with 5 μM KR for 30 min, followed by incubation with 10 μM TST for 1–6 h. (f) Cell viability MTT assay. Cells were treated as indicated. TST (10 μM) for 8 h or KR (5 μM) for 8.5 h did not influence 661W cell number. H₂O₂ (200 μM) for 6 h led to a strong decrease in cell viability, which was enhanced by pre-incubation with TST for 2 h (TST+H₂O₂). Pre-incubation with KR for 30 min followed by incubation with TST for 2 h followed by treatment with H₂O₂ for 6 h (KR+TST+H₂O₂) did not diminish the protective effect of TST. Experiments were carried out three times with similar results. Data represent the mean \pm S.D. of one representative experiment with eight replicates and are expressed as the percent of the untreated control, which was set at 100%

Image collected and cropped by CiteAb under a CC-BY license from the following publication: HDAC6 inhibition by tubastatin A is protective against oxidative stress in a photoreceptor cell line and restores visual function in a zebrafish model of inherited blindness. *Cell Death Dis* (2017)

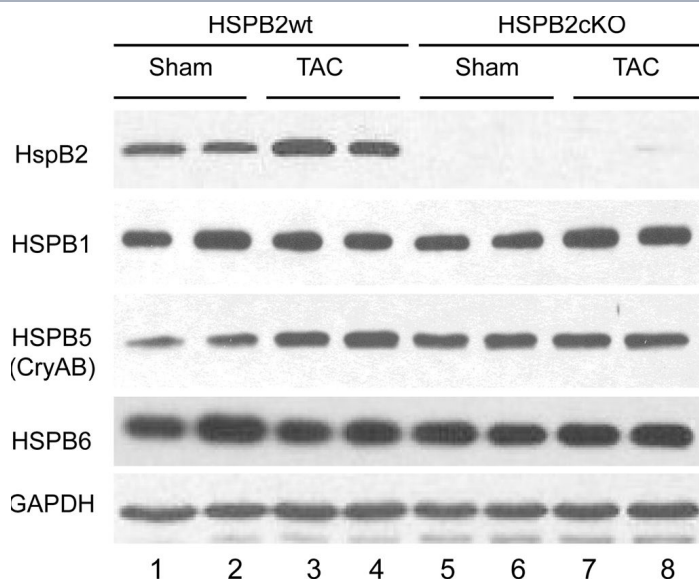


(A) Gastrocnemius muscle HSP25 and actin protein levels in Hsp25^{+/+} and Hsp25^{-/-} mice verifying the absence of HSP25. (B) HSP25 protein levels in young and old Hsp25^{+/+} mice under sedentary conditions or following 14 days of running wheel activity.

#Significantly different from sedentary ($P < 0.0001$), †significant main effect of age ($P < 0.0001$).

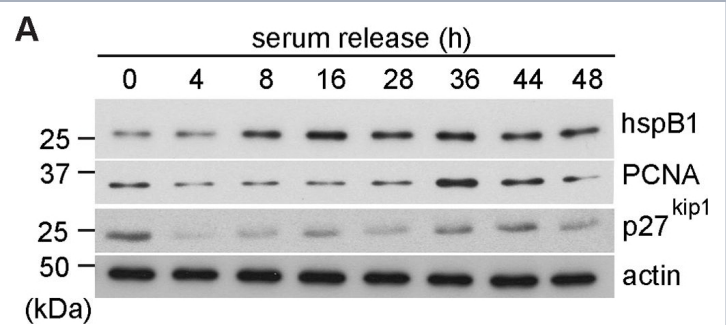
Representative western blot showing HSP25 in sedentary and running wheel young and old Hsp25^{+/+} mice. Data are presented as means \pm SE, n = 8 for all groups.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Effect of HSP25 loss on muscle contractile function and running wheel activity in young and old mice. *Front Physiol* (2014)



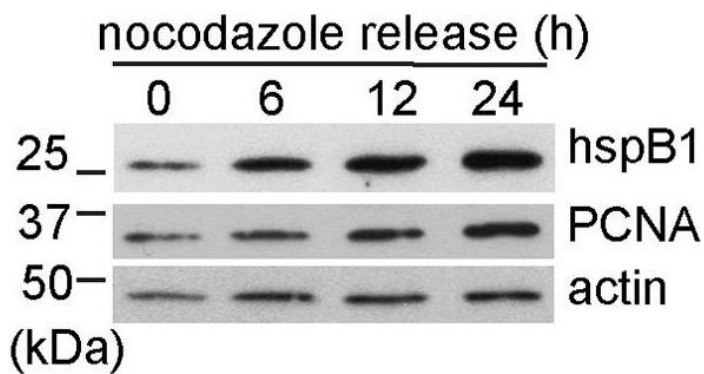
No significant changes in expression of small HSP in the absence of HSPB2 and/or under cardiac stress conditions. Hearts from HSPB2wt and HSPB2cKO were isolated at 8 weeks after sham or TAC procedure and protein extracts (soluble fraction) were analyzed by western blots probed with anti-HSPB2, HSPB1 (HSP25), HSPB5 (CRYAB) and HSPB6 (HSP20). In contrast to transcripts, protein levels were not significantly modified in samples from TAC operated mice and except for HSPB2, there was no difference between HSPB2cKO and HSPB2wt samples. Two to five samples were analyzed per group and a representative example is shown.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: HSPB2 is dispensable for the cardiac hypertrophic response but reduces mitochondrial energetics following pressure overload in mice. *PLoS One* (2012)



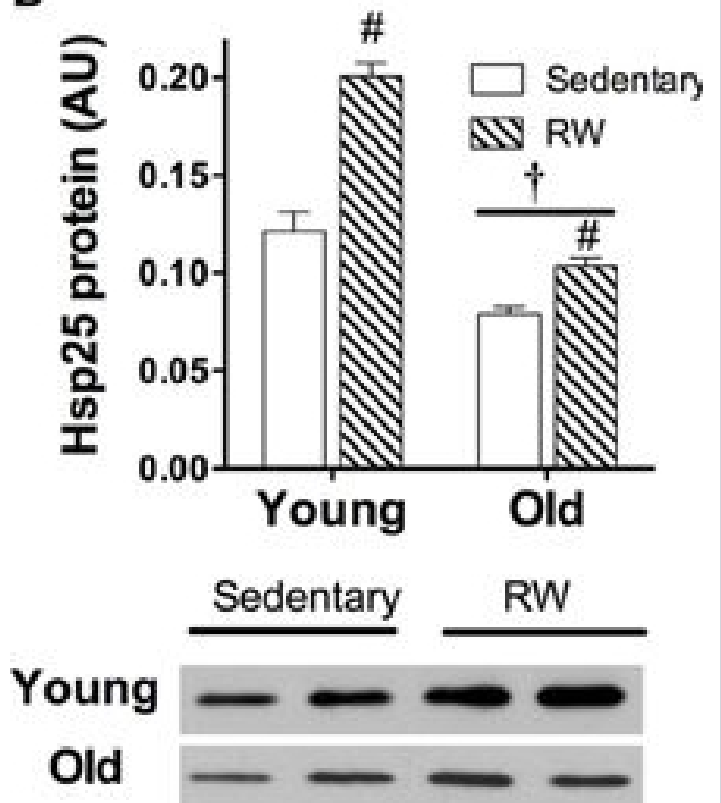
Expression of hspB1 protein and mRNA is controlled by the cell cycle. A, MEF were synchronized by serum starvation for 48h FCS-containing medium and analysed by western blot for hspB1, PCNA, p27^{kip1} and actin. B, Comparison of the expression of mRNAs for hspB1 and the cell cycle-regulated genes, Myc, and Cyclin E1 determined by qRT-PCR and normalized to GAPDH in a representative synchronized MEF serum release time course. C, Western blot for hspB1, PCNA and loading control, actin, in lysates of MEF following release from nocodazole G2/M block (40 ng/ml). All western blots shown are representative of at least two independent experiments.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Heat shock protein B1-deficient mice display impaired wound healing. *PLoS One* (2013)

C

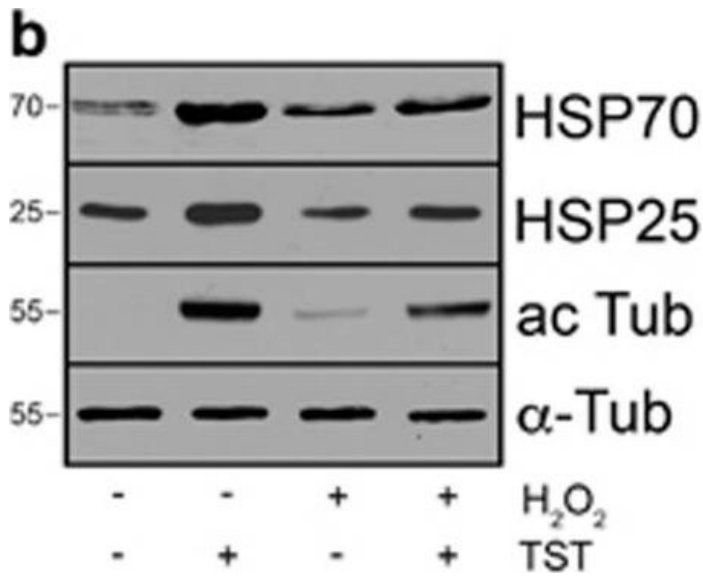
Expression of hspB1 protein and mRNA is controlled by the cell cycle. A, MEF were synchronized by serum starvation for 48h in FCS-containing medium and analysed by western blot for hspB1, PCNA, p27kip1 and actin. B, Comparison of the expression of mRNAs for hspB1 and the cell cycle-regulated genes, Myc, and Cyclin E1 determined by qRT-PCR and normalized to GAPDH in a representative synchronized MEF serum release time course. C, Western blot for hspB1, PCNA and loading control, actin, in lysates of MEF following release from nocodazole G2/M block (40 ng/ml). All western blots shown are representative of at least two independent experiments.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Heat shock protein B1-deficient mice display impaired wound healing. *PLoS One* (2013)

B

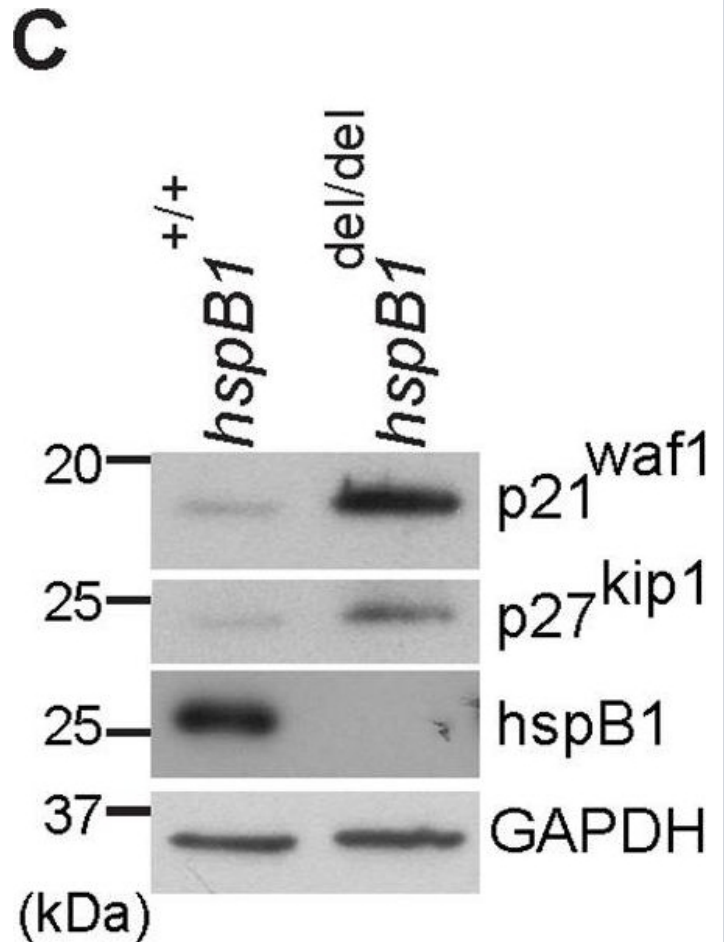
(A) Gastrocnemius muscle HSP25 and actin protein levels in Hsp25^{+/+} and Hsp25^{-/-} mice verifying the absence of HSP25. (B) HSP25 protein levels in young and old Hsp25^{+/+} mice under sedentary conditions or following 14 days of running wheel activity. #Significantly different from sedentary (P < 0.0001), †significant main effect of age (P < 0.0001). Representative western blot showing HSP25 in sedentary and running wheel young and old Hsp25^{+/+} mice. Data are presented as means ± SE, n = 8 for all groups.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Effect of HSP25 loss on muscle contractile function and running wheel activity in young and old mice. *Front Physiol* (2014)



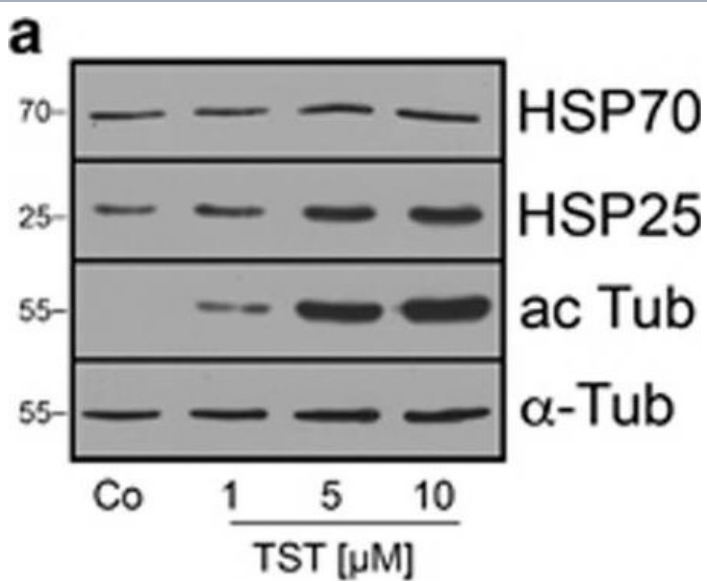
Tubastatin A induces heat-shock protein expression by activating heat-shock factor 1. 661W cells were treated with 1, 5 and 10 μ M of tubastatin A (TST) for 24 h (a), or in (b) with 10 μ M TST for 8 h, or with 200 μ M H₂O₂ for 6 h or were preincubated with 10 μ M TST for 2 h followed by incubation with 200 μ M H₂O₂ for 6 h. ac Tub, acetylated tubulin. α -Tub, α -tubulin. Co, untreated control. Quantitative evaluation of immunoblot analysis revealed a significant increase in heat-shock protein (HSP) 70 level after 8 h (c), while HSP25 was significantly enhanced after 24 h (d); n=4. (e) Heat-shock factor 1 (HSF1) activity was investigated using immunoblot analysis of 661W cell extracts that were treated 10 μ M TST for 1, 3 and 6 h, or with 5 μ M KRIBB11 (KR) for 6.5 h alone, or preincubated with 5 μ M KR for 30 min, followed by incubation with 10 μ M TST for 1–6 h. (f) Cell viability MTT assay. Cells were treated as indicated. TST (10 μ M) for 8 h or KR (5 μ M) for 8.5 h did not influence 661W cell number. H₂O₂ (200 μ M) for 6 h led to a strong decrease in cell viability, which was enhanced by pre-incubation with TST for 2 h (TST+H₂O₂). Pre-incubation with KR for 30 min followed by incubation with TST for 2 h followed by treatment with H₂O₂ for 6 h (KR+TST+H₂O₂) did not diminish the protective effect of TST. Experiments were carried out three times with similar results. Data represent the mean \pm S.D. of one representative experiment with eight replicates and are expressed as the percent of the untreated control, which was set at 100%

Image collected and cropped by CiteAb under a CC-BY license from the following publication: HDAC6 inhibition by tubastatin A is protective against oxidative stress in a photoreceptor cell line and restores visual function in a zebrafish model of inherited blindness. *Cell Death Dis* (2017)



HspB1 deficiency inhibits entry into S phase and increases the expression of p21^{waf1} and p27^{kip1}. BrdU incorporation following a 2-positive cells (mean % \pm SEM) from three independent experiments performed; **P<0.01. C, Western blot of asynchronous MEF lysates for p21^{waf1}, p27^{kip1}, hspB1 and GAPDH as a loading control with molecular weights (kDa) of markers indicated. Western blots are representative of three independent experiments.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Heat shock protein B1-deficient mice display impaired wound healing. *PLoS One* (2013)



Tubastatin A induces heat-shock protein expression by activating heat-shock factor 1. 661W cells were treated with 1, 5 and 10 μM of tubastatin A (TST) for 24 h (a), or in (b) with 10 μM TST for 8 h, or with 200 μM H₂O₂ for 6 h or were preincubated with 10 μM TST for 2 h followed by incubation with 200 μM H₂O₂ for 6 h. ac Tub, acetylated tubulin. α -Tub, α -tubulin. Co, untreated control. Quantitative evaluation of immunoblot analysis revealed a significant increase in heat-shock protein (HSP) 70 level after 8 h (c), while HSP25 was significantly enhanced after 24 h (d); n=4. (e) Heat-shock factor 1 (HSF1) activity was investigated using immunoblot analysis of 661W cell extracts that were treated 10 μM TST for 1, 3 and 6 h, or with 5 μM KRIBB11 (KR) for 6.5 h alone, or preincubated with 5 μM KR for 30 min, followed by incubation with 10 μM TST for 1–6 h. (f) Cell viability MTT assay. Cells were treated as indicated. TST (10 μM) for 8 h or KR (5 μM) for 8.5 h did not influence 661W cell number. H₂O₂ (200 μM) for 6 h led to a strong decrease in cell viability, which was enhanced by pre-incubation with TST for 2 h (TST+H₂O₂). Pre-incubation with KR for 30 min followed by incubation with TST for 2 h followed by treatment with H₂O₂ for 6 h (KR+TST+H₂O₂) did not diminish the protective effect of TST. Experiments were carried out three times with similar results. Data represent the mean \pm S.D. of one representative experiment with eight replicates and are expressed as the percent of the untreated control, which was set at 100%

Image collected and cropped by CiteAb under a CC-BY license from the following publication: HDAC6 inhibition by tubastatin A is protective against oxidative stress in a photoreceptor cell line and restores visual function in a zebrafish model of inherited blindness. *Cell Death Dis* (2017)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name HspB1, Heat shock protein 25

Application Electron microscopy, ICC, IF, IP, WB

Application Notes Detects a band of ~25kDa by Western blot.

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

GenBank ID L07577

Host Rabbit

Immunogen Recombinant mouse Hsp25.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Immunoprecipitation (1:100)Western Blot (1:1,000, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source Purified from rabbit serum.

Species Reactivity Bovine, Dog, Gerbil, Guinea pig, Hamster, Human, Mouse, Porcine, Rat

UniProt ID P14602

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



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)