

# HSF1 monoclonal antibody (10H8)

HSFs (Heat Shock family of transcription factors), which consists of HSF 1-4, bind to highly conserved Heat shock elements (HSEs) in the promoter regions of heat shock genes, ultimately regulating the expression of Heat shock proteins (Hsps). On exposure to heat shock and other stresses, HSF1 localizes within seconds to discrete nuclear granules and on recovery from stress, HSF1 rapidly dissipates from the stress granules to a diffuse nucleoplasmic distribution.

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

Citations: 12

[View Online »](#)

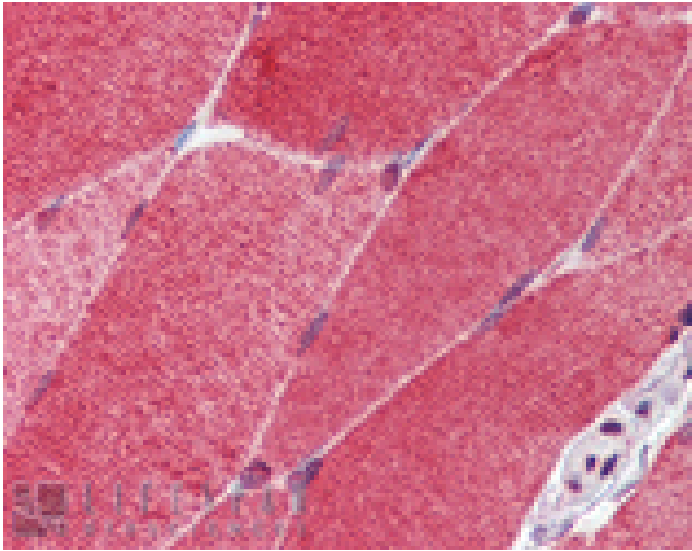
## Ordering Information

[Order Online »](#)

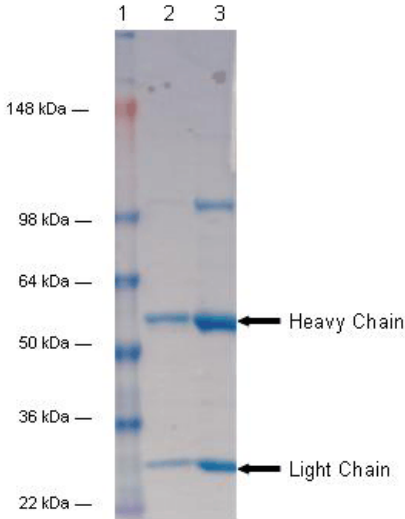
ADI-SPA-950-D	50µg
ADI-SPA-950-F	200µg

## Manuals, SDS & CofA

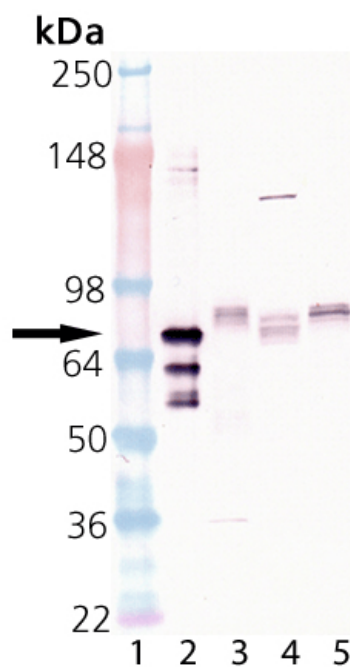
[View Online »](#)



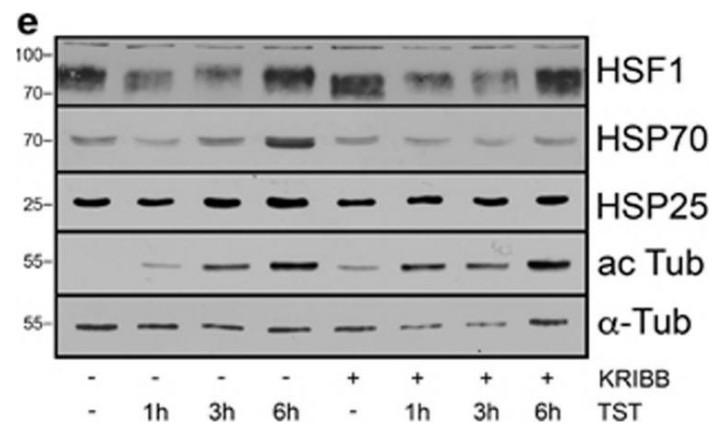
Immunohistochemistry analysis of human skeletal muscle tissue stained with HSF1, mAb (10H8) at 10µg/ml.



SDS-PAGE analysis of ADI-SPA-950: Lane 1: MW Marker, Lane 2: 1µg and Lane 3: 5µg of HSF1 monoclonal antibody (10H8), resolved under partially reduced condition.



Western Blot analysis of ADI-SPA-950: Lane 1: MW Marker, Lane 2: HSF1 (human), (recombinant) (His-tag), (Prod. No. ADI-SPP-900), Lane 3: HeLa lysate, (Prod. No. ADI-LYC-HL100), Lane 4: HeLa (heat shocked) lysate, (Prod. No. ADI-LYC-HL102), Lane 5: PC-12 lysate, (Prod. No. ADI-LYC-PC100).



Tubastatin A induces heat-shock protein expression by activating heat-shock factor 1. 661W cells were treated with 1, 5 and 10  $\mu\text{M}$  of tubastatin A (TST) for 24 h (a), or in (b) with 10  $\mu\text{M}$  TST for 8 h, or with 200  $\mu\text{M}$  H<sub>2</sub>O<sub>2</sub> for 6 h or were preincubated with 10  $\mu\text{M}$  TST for 2 h followed by incubation with 200  $\mu\text{M}$  H<sub>2</sub>O<sub>2</sub> for 6 h. ac Tub, acetylated tubulin.  $\alpha$ -Tub,  $\alpha$ -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  $\mu\text{M}$  TST for 1, 3 and 6 h, or with 5  $\mu\text{M}$  KRIBB11 (KR) for 6.5 h alone, or preincubated with 5  $\mu\text{M}$  KR for 30 min, followed by incubation with 10  $\mu\text{M}$  TST for 1–6 h. (f) Cell viability MTT assay. Cells were treated as indicated. TST (10  $\mu\text{M}$ ) for 8 h or KR (5  $\mu\text{M}$ ) for 8.5 h did not influence 661W cell number. H<sub>2</sub>O<sub>2</sub> (200  $\mu\text{M}$ ) for 6 h led to a strong decrease in cell viability, which was enhanced by pre-incubation with TST for 2 h (TST+H<sub>2</sub>O<sub>2</sub>). Pre-incubation with KR for 30 min followed by incubation with TST for 2 h followed by treatment with H<sub>2</sub>O<sub>2</sub> for 6 h (KR+TST+H<sub>2</sub>O<sub>2</sub>) 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** HSTF1, Heat shock factor protein 1

**Application** IHC (PS), WB

**Application Notes** Predicted MW of ~60kDa. Detects a band of ~85kDa by Western blot.

**Clone** 10H8

**Formulation** Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

**Gene/Protein Identifier** NM\_008296 (RefSeq)

**Host** Rat

**Immunogen** Recombinant mouse HSF1.

**Isotype** IgG1

**Purity Detail** Protein G affinity purified.

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

**Source** Purified from ascites.

**Species Reactivity** Human, Monkey, Mouse, Rabbit, Rat



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)

**Worry-free Guarantee**

**This antibody is covered by our Worry-Free Guarantee.**

For local distributors and detailed product information visit us online: [enzo.com](http://enzo.com) All product names, logos, and brands are the property of their respective owners.