# 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 »

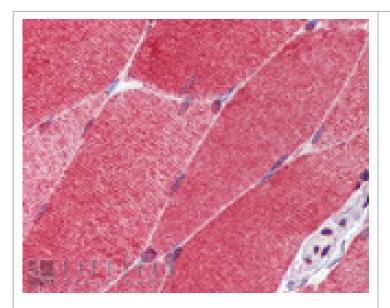
# **Ordering Information**

Order Online »

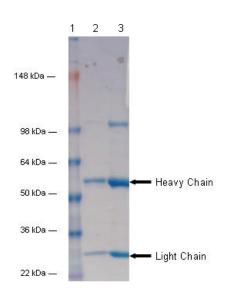
ADI-SPA-950-D	50µg
ADI-SPA-950-F	200μg

Manuals, SDS & CofA

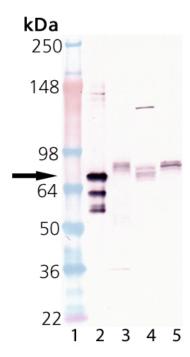
**View Online** »



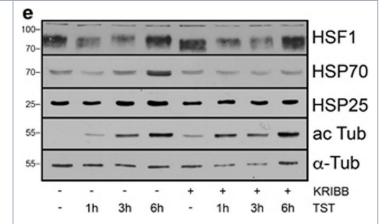
Immunohistochemistry analysis of human skeletal muscle tissue stained with HSF1, mAb (10H8) at  $10\mu g/ml$ .



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 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.  $\alpha$ -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) Heatshock 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  $\mu$ M KR for 30 min, followed by incubation with 10  $\mu$ M 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±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.

lsotype lgG1

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions 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 ascites.

Species Reactivity Human, Monkey, Mouse, Rabbit, Rat

**Worry-free Guarantee** 

This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

