

# HSP20 polyclonal antibody

Hsp20 is a member of the family of small heat shock proteins (sHsps), with homology to Hsp25, Hsp27, and alphaB-crystallin. Highest expression of Hsp20 is found in skeletal, smooth, and cardiac muscle. Under heat shock, Hsp20 translocates from the cytosol to the nucleus, although the amount of Hsp20 in the cell does not increase like other Hsps. Hsp20 oligomerization and interaction with polymeric actin filaments is regulated via phosphorylation by PKA and PKG.

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

Citations: 2

[View Online »](#)

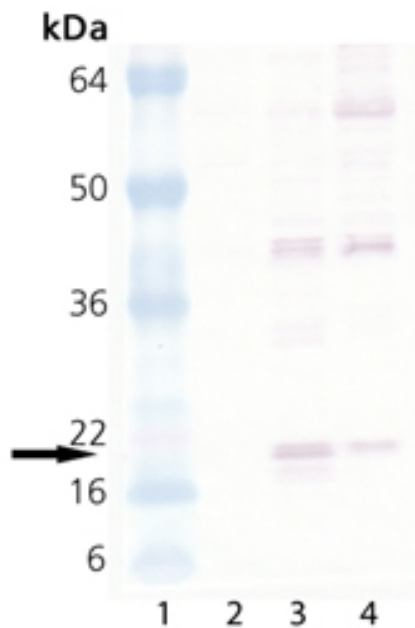
Ordering Information

[Order Online »](#)

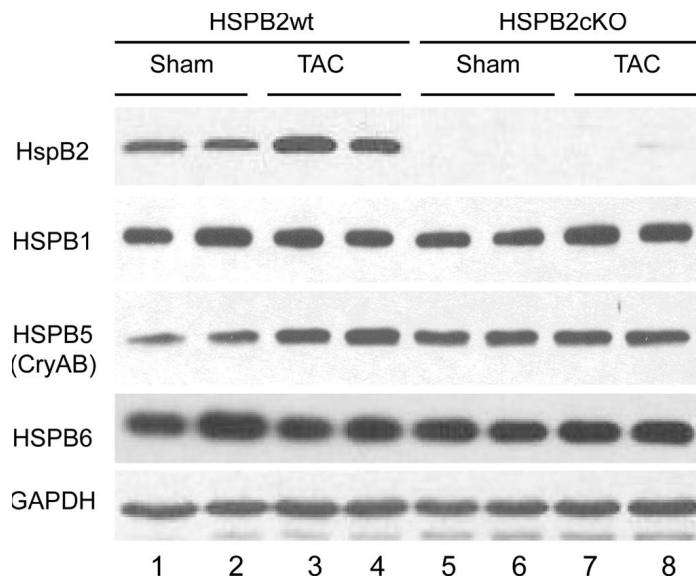
ADI-SPA-796-F	200µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of HSP20, pAb (Prod. no. ADI-SPA-796): Lane 1: MW Marker, Lane 2: HeLa (heat shocked) (Prod. No. ADI-LYC-HL101), Lane 3: 3T3 (heat shocked) (Prod. No. ADI-LYC-3T101), Lane 4: PC-12 (heat shocked) (Prod. No. ADI-LYC-PC101)



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)

## 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	HspB6, Heat shock protein 20
Application	WB
Application Notes	Detects a band of ~20kDa by Western blot.
Formulation	Liquid. In 0.1M phosphate buffer, pH 7.0, containing 3% BSA and 0.09% sodium azide.
Host	Rabbit
Immunogen	Rat skeletal muscle Hsp20.
Purity Detail	Peptide 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 rabbit serum.
Species Reactivity	Bovine, Dog, Fish, Human, Mouse, Rat
UniProt ID	P97541

**Worry-free Guarantee**

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



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-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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

Last modified: May 29, 2024