

[pSer326]HSF1 polyclonal antibody

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 »](#)

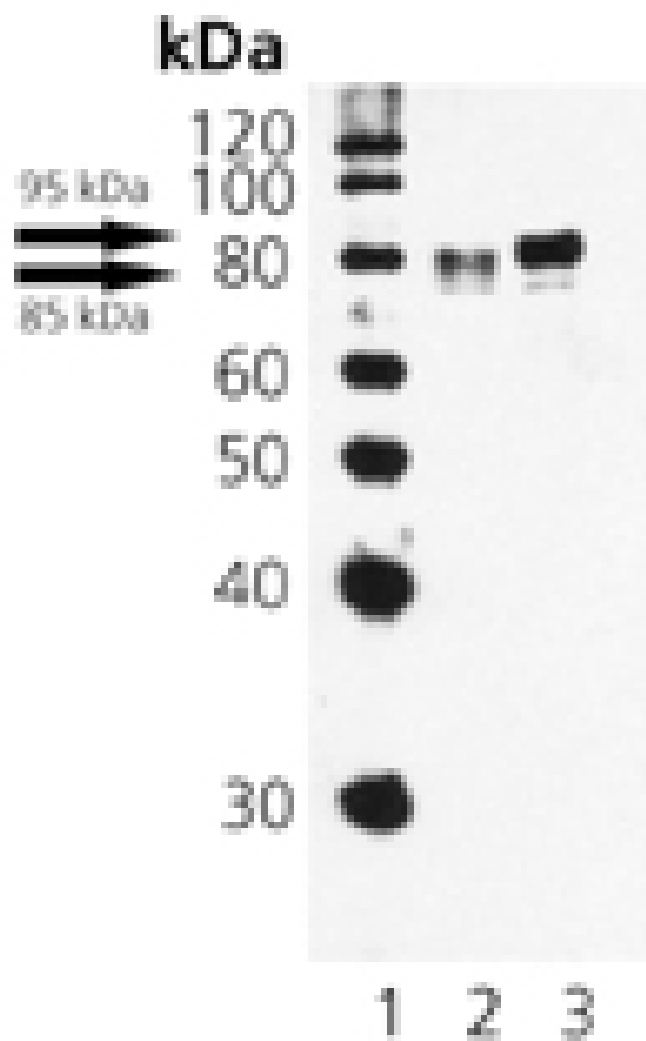
Ordering Information

[Order Online »](#)

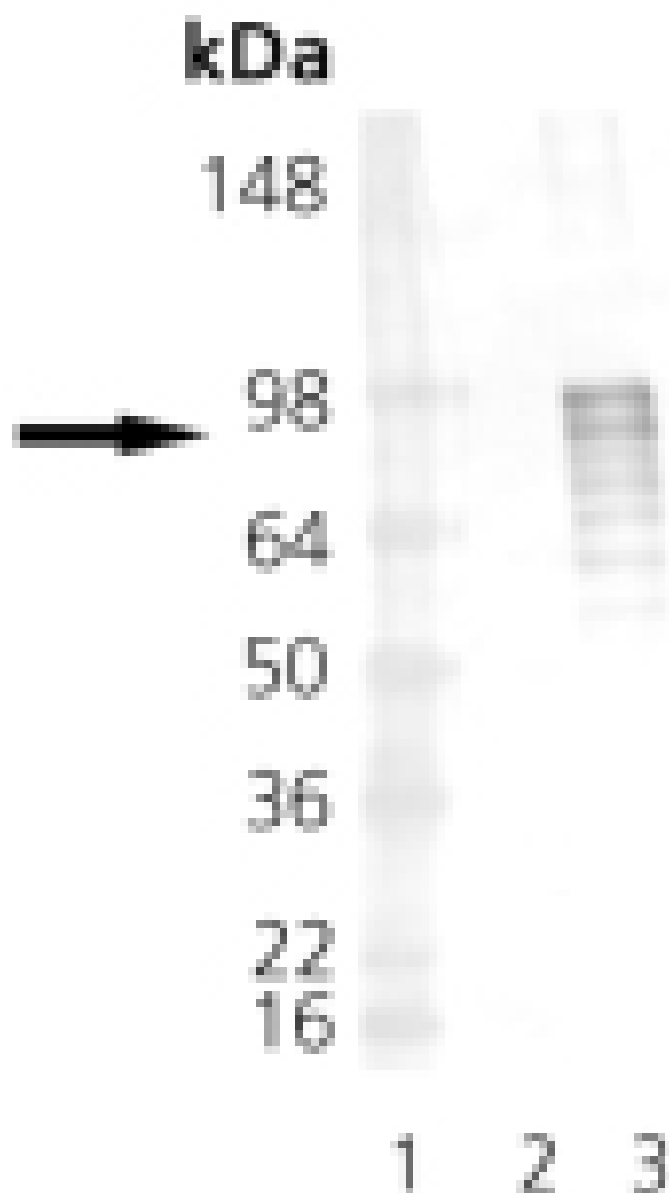
ADI-SPA-902-D	50µg
ADI-SPA-902-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of HSF1 (pSer326): Lane 1: MW marker, Lane 2: HeLa, Lane 3: HeLa, Heat-Shocked (No recovery time), probed with HSF1 (pSer326) pAb.



Western Blot analysis of Recombinant HSF1 and HSF1 (phospho): Lane 1: MWM, Lane 2: Recombinant HSF1, Lane 3: Recombinant HSF1 (phospho), probed with HSF1 (pSer326) pAb.

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Alternative Name	HSTF1, Heat shock factor protein 1
Application	WB
Application Notes	Predicted MW of ~60kDa. Detects a band of 85-95kDa by Western blot depending on hyperphosphorylation of the native protein.
Formulation	Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to the N-terminus of human HSF1 phosphorylated at Ser 326.
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:1,000, ECL)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	Human
UniProt ID	Q00613
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-
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)