

[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.

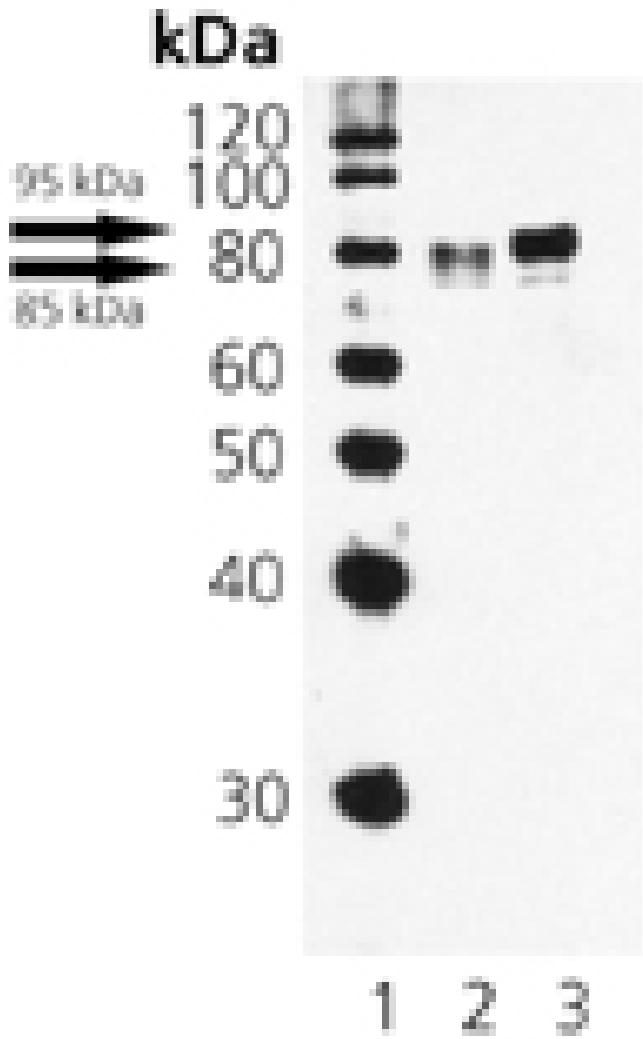
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 12 [View Online »](#)

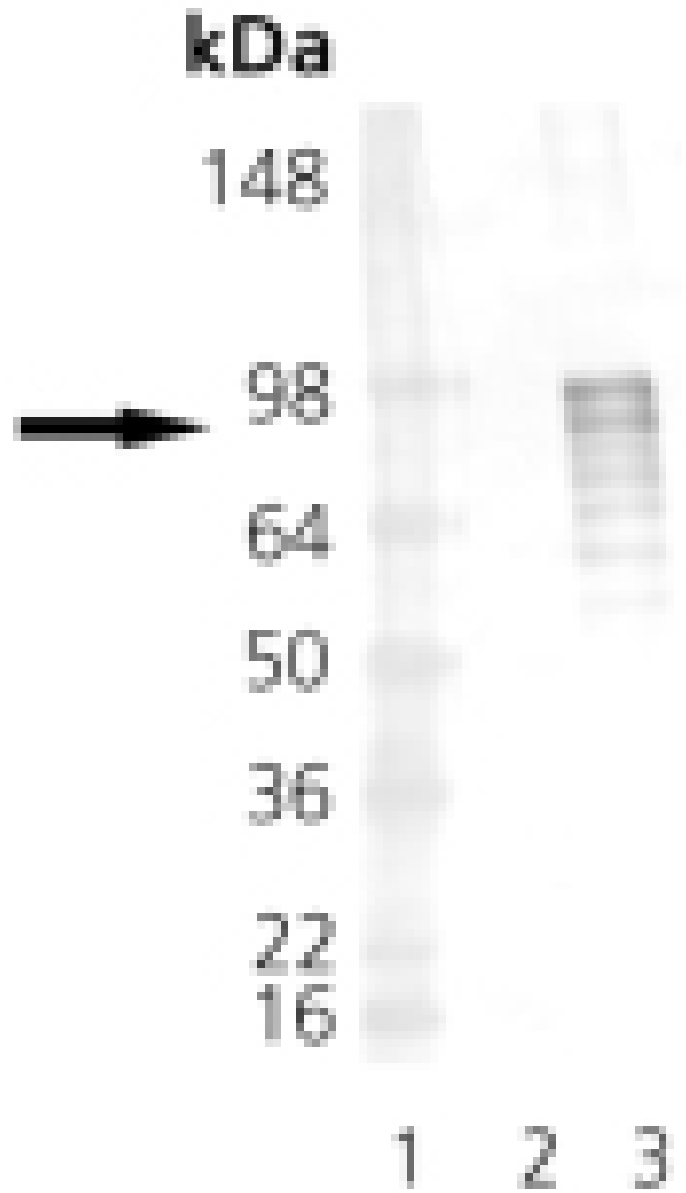
Ordering Information [Order Online »](#)

ADI-SPA-902-D	50µg
---------------	------

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 & 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 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](#)

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)