

HSF1 (human), (recombinant) (His- tag)

In response to adverse changes in their environment, cells from many organisms increase the expression of a class of proteins referred to as heat shock or stress proteins. The expression of heat shock proteins is mediated by the binding of a heat shock transcriptional factor (HSF) to a heat shock response element (HSE) which is located upstream of heat shock protein genes. Two heat shock transcriptional factors, termed HSF1 and HSF2, have been identified in mice and humans. In unstressed cells, HSF1 is a non-DNA binding monomeric protein present in the cytosol and nuclear compartments. However, in response to heat and heavy metals, HSF1 rapidly converts to a transcriptionally active form - a process requiring oligomerization to a trimer, serine phosphorylation, and translocation to the nucleus. HSF2, on the other hand, is not activated by heat, but is involved in regulating heat shock gene transcription during cell differentiation and development.

Citations: 4

[View Online »](#)

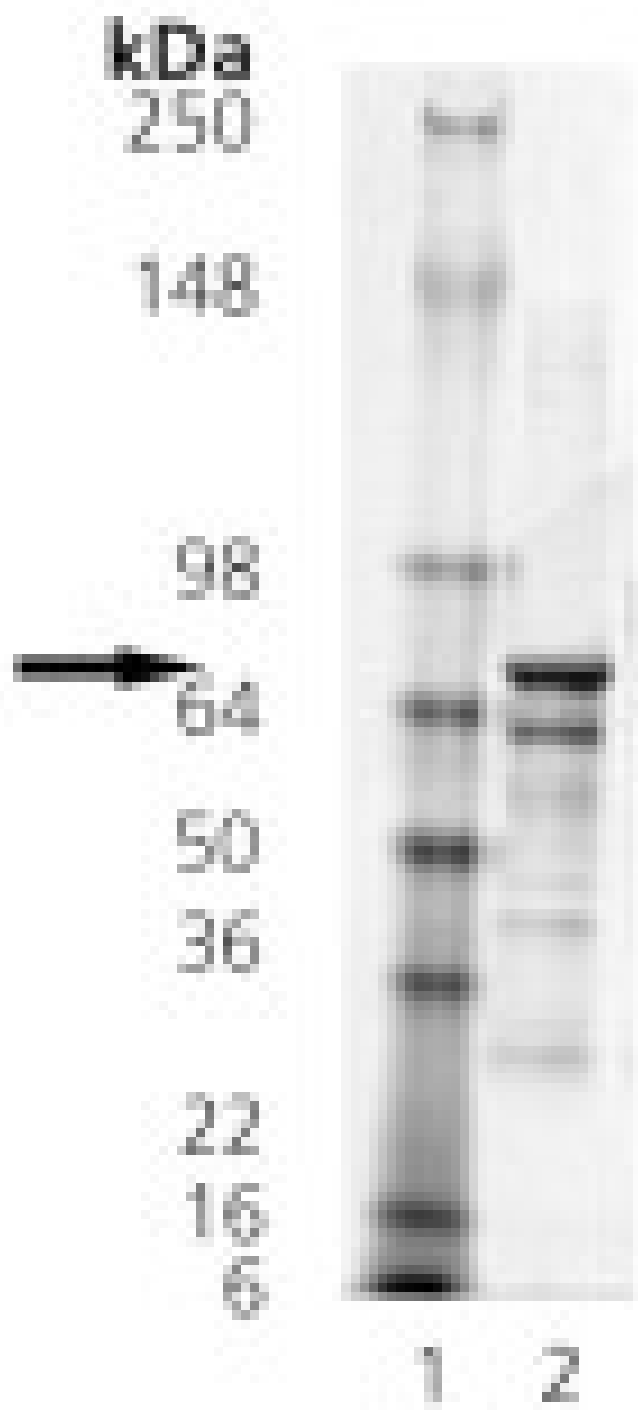
Ordering Information

[Order Online »](#)

ADI-SPP-900-F	200µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis: Lane 1: MWM, Lane 2: 2 μ g HSF1.



Western Blot analysis: Lane 1: MW marker, Lane 2: HSF1 probed with HSF1 pAb.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	HSTF1, Heat shock factor protein 1
Application Notes	Western blot control.
Formulation	Liquid. In 1x TRIS buffered saline containing 5% glycerol.
MW	~60kDa (predicted), ~70kDa (observed due to His-tag)
Purity	≥90% (~60% full-length)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> . Human HSF1 is fused at the N-terminus to a His-tag.
UniProt ID	Q00613

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)