

HSF1 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog # H25-30G

Lot # F385-4

Product Description

Recombinant full length human HSF1 was expressed in E.coli cells using an N-terminal GST tag. The gene accession number is [NM_005526](#).

Gene Aliases

HSTF1

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

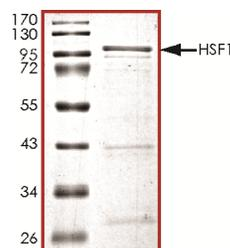
Scientific Background

HSF1 is a member of the heat shock transcription factor family. Protein-damaging stress lead to the activation of HSF1 which binds to upstream regulatory sequences in the promoters of heat shock genes leading to enhanced heat shock gene expression (1). The activation of HSF1 proceeds through a multi-step pathway, involving a monomer-to-trimer transition, nuclear accumulation and extensive posttranslational modifications. HSF1 activity is regulated at different levels by heat shock proteins and co-chaperones and is modulated further by a number of mechanisms involving other stress-regulated aspects of cell metabolism (2).

References

1. Pirkkala, L. et al: Roles of the heat shock transcription factors in regulation of the heat shock response and beyond. *FASEB J.* 2001 May;15(7):1118-31.
2. Voellmy, R.: Feedback regulation of the heat shock response. *Handb Exp Pharmacol.* 2006;(172):43-68.

Purity



The purity of HSF1 was determined to be **>85%** by densitometry. Approx. MW **96kDa**.

Related Protein Products

HSP10	Cat # H29-54G
HSP27	Cat # H30-54G
HSP40	Cat # H32-54G
HSP70	Cat # H34-54G
HSP70	Cat # H34-54H
HSP90β	Cat # H36-54H

HSF1 Protein

Full-length recombinant protein expressed in E.coli cells

Catalog Number	H25-30G
Specific Lot Number	F385-4
Purity	>85%
Concentration	0.2µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.