

Phospho-HSF1 (S326) Recombinant Monoclonal Antibody [2A7]

(Catalog # A74142)

Background

Functions as a stress-responsive transcription factor, orchestrating the heat shock response (HSR) by activating heat shock proteins (HSPs) that safeguard cells from stress-induced damage. In unstressed cells, resides in an HSP90-containing complex, remaining inactive. Upon stress, it trimerizes, binds specific heat shock elements (HSEs) in HSP gene promoters, and activates their transcription. This process is reversible. Binds chromatin at heat shock gene promoters and has non-transcriptional roles too. Represses Ras-induced c-fos transcription during heat stress, aids HSP70 mRNA processing and export, regulates mitotic progression, and negatively influences DNA repair. Facilitates stress-driven cancer cell proliferation through IER5 dependence.

Description

Phospho-HSF1 (S326) Recombinant Monoclonal Antibody [2A7]. Unconjugated. Raised in: HEK293F Cell.

Formulation

Buffer: Rabbit IgG in phosphate buffered saline, pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.

Specificity

Human

Isotype

Rabbit IgG

Uniprot ID

Q00613

Purification

Affinity Chromatography

Immunogen

A synthesized peptide derived from human Phospho-HSF1 (S326)

Storage

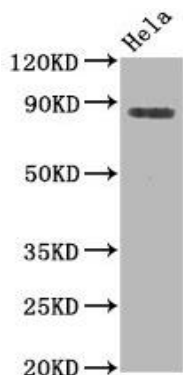
Upon receipt, store at -20°C or -80°C. Avoid repeated freeze.

Alternative Names

Heat shock factor protein 1, HSF 1, Heat shock transcription factor 1, HSF1

Application

ELISA, WB, IHC; Recommended dilution: WB:1:500-1:5000, IHC:1:50-1:200



Western Blot

Positive WB detected in HeLa whole cell lysate

All lanes Phospho-HSF1 antibody at 0.8ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 82 KDa

Observed band size: 82 KDa