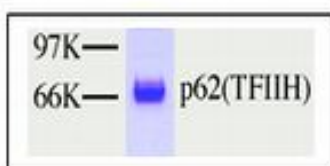


TFIIH-p62

Transcription Factor IIH, p62 subunit
human, recombinant, *E. coli*

Cat. No.	Amount
PR-711	10 µg



For *in vitro* use only
Quality guaranteed for 12 months
Store at -80°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 20 mM Tris-HCl pH 8.0, 100 mM KCl, 0.2 mM EDTA, 1 mM DTT, 20% glycerol.

Activity

100 ng are sufficient for a protein-protein interaction assay.

Molecular Weight

68 kDa

Purity

> 95% by SDS-PAGE

Description

TFIIH, a multisubunit complex is involved in several biological fundamental mechanisms of the cell: transcription, nucleotide excision repair, and cell cycle regulation. p62 is one of the six subunits that constitutes the core of TFIIH. Analysis of the expression of the p62 gene reveals an over-expression in testis tissue. This subunit of TFIIH participates in a variety of protein-protein interactions. For example, Rb competes with TBP and p62 for binding to E2F thus repressing E2F-mediated transactivation, herpes simplex virus VP16 and human p53, directly interact with the p62 subunit of TFIIH. In addition, TFIIH, via p62 phosphorylation is the major target for mitotic inactivation of transcription.

Recombinant p62 protein is isolated from an *E. coli* strain that carries the coding sequence of human p62 under the control of a T7 promoter.

p62 has been used for protein-protein interactions assays.

Protein is greater than 95% homogeneous and contains no detectable protease, DNase, and RNase activity.

Selected References:

- Feaver *et al.* (1991) Purification and characterization of yeast RNA polymerase II transcription factor b. *J. Biol. Chem.* **266**:19000.
- Flores *et al.* (1992) Factors involved in specific transcription by mammalian RNA polymerase II. Identification and characterization of factor IIH. *J. Biol. Chem.* **267**:2786.
- Fischer *et al.* (1992) Cloning of the 62-kilodalton component of basic transcription factor BTF2. *Science* **257**:1392.
- Perez *et al.* (1998) Genomic organization and promoter characterization of the mouse and human genes encoding p62 subunit of the transcription/DNA repair factor TFIIH. *Gene* **213**:73.
- Pearson *et al.* (1997) Modular organization of the E2F1 activation domain and its interaction with general transcription factors TBP and TFIIH. *Oncogene* **15**:2643.
- Xiao *et al.* (1994) Binding of basal transcription factor TFIIH to the acidic activation domains of VP16 and p53. *Mol. Cell. Biol.* **14**:7013.
- Long *et al.* (1998) Repression of TFIIH transcriptional activity and TFIIH-associated cdk7 kinase activity at mitosis. *Mol. Cell. Biol.* **18**:1467.