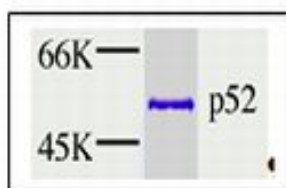


p52

Transcriptional Coactivator
human, recombinant, *E. coli*

Cat. No.	Amount
PR-728	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 and 20% glycerol.

Activity

1 ng is sufficient for a gel mobility shift assay in a 20 µl reaction, 20 ng are sufficient for reconstituted transcription assay and 100 ng are sufficient for a protein-protein interaction assay.

Molecular Weight

54 kDa

Purity

> 95% by SDS-PAGE

Description

The human p52 protein is a non-TAF transcription coactivator that mediates activator-dependent transcription by RNA Polymerase II. The function of p52 is through interactions with transcriptional activators and the basal transcription machinery. In addition, p52 could also interact with several cellular proteins including the transcription coactivator PC4, the essential splicing factor ASF/SF2 and the nuclear protein nucleolin.

Recombinant p52 protein (wild type, 333 amino acids) is isolated from an *E. coli* strain that carries the coding sequence of human p52 under the control of T7 promoter and purified by the combination of an affinity column and FPLC chromatography.

Recombinant p52 has been utilized for *in vitro* function studies, including transcription, splicing, protein-DNA/RNA and protein-protein interactions.

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

Selected References:

Ge *et al.* (1998) Isolation of cDNAs encoding novel transcription coactivators p52 and p75 reveals an alternate regulatory mechanism of transcriptional activation. *EMBO J.* **17**:6723.

Ge *et al.* (1998) A novel transcriptional coactivator, p52, functionally interacts with the essential splicing factor ASF/SF2. *Molecular Cell* **2**:751.

Ge (2000) UPA, a universal protein array system for quantitative detection of protein-protein, protein-DNA, protein-RNA and proteinligand interactions. *Nucleic Acid Res.* **28**:e3.