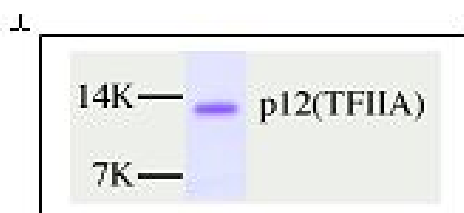


## TFIIA-p12

Transcription Factor IIA, p12 subunit  
human, recombinant, *E. coli*

Cat. No.	Amount
PR-701	10 $\mu$ g



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 of p12 is sufficient for a gel mobility shift assay, 10 ng are sufficient for transcription assay and 100 ng are sufficient for a protein-protein interaction assay.

### Application

Recombinant p12 in combination with other subunits of TFIIA has been applied for *in vitro* transcription assays, gel mobility shift assays and protein-protein interaction assays.

### Molecular Weight

12 kDa

### Purity

> 95% by SDS-PAGE

### Description

p12 is a small subunit of the transcription factor IIA and has been shown to be required for both basal and activated transcription.

Recombinant p12, along with two other subunits ( $\alpha$  and  $\beta$ ) can potentiate transcriptional activation, whereas p12 along with  $\beta$ -subunit is able to function in an antirepression.

p12 subunit of TFIIA is isolated from a strain of *E. coli* that contains the coding sequence for human TFIIA p12 under the control of T7 promoter.

### Selected References:

- DeJong *et al.* (1995) Human general transcription factor TFIIA: characterization of a cDNA encoding the small subunit and requirement for basal and activated transcription. *Proc. Natl. Acad. Sci. USA* **92**:3313.
- Ozer *et al.* (1994) Molecular cloning of the small ( $\gamma$ ) subunit of human TFIIA reveals functions critical for activated transcription. *Genes & Dev.* **8**:2324.
- Sun *et al.* (1994) Reconstitution of human TFIIA activity from recombinant polypeptides: a role in TFIID-mediated transcription. *Genes & Dev.* **8**:2336.
- Ma *et al.* (1996) Separation of the Transcriptional Coactivator and Antirepression Functions of Transcription Factor IIA. *Proc. Natl. Acad. Sci. USA* **93**:6583.