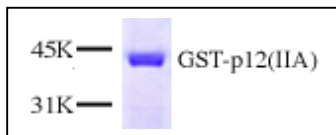


TFIIA^{GST} (p12)

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

Cat. No.	Amount
PR-786	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

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

Purity

> 95% by SDS-PAGE

Description

p12 is a small subunit (γ) of the Transcription Factor IIA and has been shown to be required for both basal and activated transcription. Liquid. Recombinant p12, along with two other subunits (α and β) can potentiate transcriptional activation, whereas p12 along with β -subunit is able to function in an antirepression. GST-p12 is isolated from a strain of *E. coli* that contains the coding sequence for human TFIIA γ -subunit under the control of a 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.
- Kaludov *et al.* (2000) MeCP2 driven transcriptional repression *in vitro*: selectivity for methylated DNA, action at a distance and contacts with the basal transcription machinery. *Nucleic Acids Res.* **28**:1921.