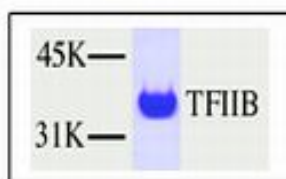


TFIIB

Transcription Factor IIB

human, recombinant, *E. coli*

Cat. No.	Amount
PR-702	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

1 ng is sufficient for gel mobility shift assay in a 20 µl reaction to super-shift TBP-DNA complex, 10 ng are sufficient for a 25 µl reconstituted transcription assay and 100 ng are sufficient for a protein-protein interaction assay detected by immuno-blot system.

Molecular Weight

36 kDa

Purity

> 95% by SDS-PAGE

Description

The transcription factor IIB (TFIIB) is an essential factor for transcription by RNA Polymerase II. TFIIB has been shown to be required for selective binding of RNA Polymerase II, TFIIF and TFIID-DNA complex, and for specifying the start site of transcription. Human TFIIB is a single, 36 kDa polypeptide, homologous to yeast factor ϵ , the product of the SUA7 gene.

Isolated from an *E. coli* strain that carries the coding sequence for human TFIIB under the control of T7 promoter. Recombinant TFIIB has been used for *in vitro* transcription assays, gel mobility shift assay.

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

Selected References:

- Malik *et al.* (1991) Sequence of general transcription factor TFIIB and relationships to other initiation factors. *Proc. Natl. Acad. Sci. USA* **88**:9553.
Conaway *et al.* (1993) General initiation factors for RNA polymerase II. *Annu. Rev. Biochem.* **62**:161.
Orphanides *et al.* (1996) The general transcription factors of RNA polymerase II. *Genes & Dev.* **10**:2657.
Buratowski *et al.* (1989) Five intermediate complexes in transcription initiation by RNA polymerase II. *Cell* **56**:549.
Ge *et al.* (1996) *Methods Enzymol.* **274**:57.