

## TFIIB<sup>GST</sup>

### Transcription Factor IIB

human, recombinant, *E. coli*

Cat. No.	Amount
PR-787	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 detected by immuno-blot system.

#### 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, 35 kDa polypeptide, homologous to yeast factor  $\epsilon$ , the product of the SUA7 gene.

The Transcription Factor IIB (TFIIB) is an essential factor for transcription by RNA polymerase II.

Isolated from an *E. coli* strain that carries the coding sequence for human TFIIB under the control of a T7 promoter.

GST-TFIIB has been used for protein-protein interaction assays.

#### 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.
- 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.