

TBP^{GST}

TATA box Binding Protein

human, recombinant, *E. coli*

Cat. No.	Amount
PR-781	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

100 ng of the protein is sufficient for a protein-protein interaction assay detected by immuno-blot system.

Purity

>95% by SDS-PAGE

Description

The TATA-binding Protein (TBP) is believed to function as an essential factor of the general transcription machinery and to be involved in transcription by all three eukaryotic RNA Polymerases (Pol I, II, and III). TBP specifically binds to TATA element at the promoter region and interacts with numerous transcription factors, including TBP-associated Factors (TAFs), activators, and some tumor suppressor proteins.

Isolated from an *E. coli* strain that carries the coding sequence for human TBP under the control of a T7 promoter. GST-TBP can be used for protein-protein interaction assay.

Selected References:

- Horikoshi *et al.* (1988) Transcription factor ATF interacts with the TATA factor to facilitate establishment of a preinitiation complex. *Cell* **54**:1033.
- Hernandez *et al.* (1993) TBP, a universal eukaryotic transcription factor? *Genes & Dev.* **7**:1291.
- 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.