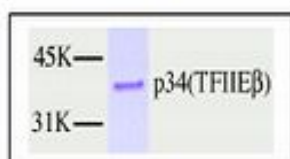


TFII β , p34

Transcription Factor IIE, β -subunit
human, recombinant, *E. coli*

Cat. No.	Amount
PR-705	10 μ 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

20 ng are sufficient for an *in vitro* reconstituted transcription assay in the presence of 56 kDa subunit, 100 ng are sufficient for a protein-protein interaction assay.

Molecular Weight

36 kDa

Purity

> 95% by SDS-PAGE

Description

The human Transcription Factor IIE (TFII β) is composed of 56 kDa (α) and 34 kDa (β) subunits and is shown to be a heterotetramer.

TFII β binds to RNA Polymerase II in solution and joins the preinitiation complex probably concomitant with RNA Polymerase II and TFII α .

The recombinant protein is purified from an *E. coli* strain that contains the coding sequence of human TFII β subunit under the control of T7 promoter.

Both subunits are required to reconstitute the functional transcription factor IIE.

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

Selected References:

Ohkuma *et al.* (1990) Factors involved in specific transcription by mammalian RNA polymerase II: purification and characterization of general transcription factor TFII β . *Proc. Natl. Acad. Sci. USA* **87**:9163.

Inostroza *et al.* (1991) Factors involved in specific transcription by mammalian RNA polymerase II. Purification and functional analysis of general transcription factor IIE. *J. Biol. Chem.* **266**:9304.

Flores *et al.* (1990) Factors involved in specific transcription by mammalian RNA polymerase II. Purification and subunit composition of transcription factor IIF. *J. Biol. Chem.* **265**:5629.

Maldonado *et al.* (1996) Purification of human RNA polymerase II and general transcription factors. *Methods Enzymol.* **274**:72.