

TFIIA^{GST} (p19)

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

Cat. No.	Amount
PR-785	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

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

Purity

> 95% by SDS-PAGE

Description

The Transcription Factor IIA (TFIIA) has been shown to bind to the TBP-DNA complex and to increase the affinity of TBP for the TATA element. Human TFIIA consists of three subunits of 35 kDa (α -subunit), 19 kDa (β -subunit) and 12 kDa (γ -subunit). The α - and β - subunits are derived from the product, p55, of a single gene by an unknown mechanism. However, recombinant p55, in combination with a 12 kDa subunit (γ -subunit), retains native TFIIA activity.

GST-TFIIA(p19) is isolated from an *E. coli* strain that carries the coding sequence of TFIIA γ -subunit under the control of a T7 promoter.

GST-TFIIA (p19) has been applied in protein-protein interaction assays.

Selected References:

- Buratowski *et al.* (1989) Five intermediate complexes in transcription initiation by RNA polymerase II. *Cell* **56**:549.
Ranish *et al.* (1991) The yeast general transcription factor TFIIA is composed of two polypeptide subunits. *J. Biol. Chem.* **266**:19320.
DeJong *et al.* (1993) A single cDNA, hTFIIA/alpha, encodes both the p35 and p19 subunits of human TFIIA. *Genes & Dev.* **7**:2220.