

α -TFIIIE β (anti-p34)

anti-p34 subunit of human Transcription Factor II E β

rabbit, polyclonal

Cat. No.	Amount
ABD-004	50 μ g

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied in PBS.

Description

The human Transcription Factor IIE (TFIIE) is composed of 56 kDa (α) and 34 kDa (β) subunits and is shown to be a heterotetramer. The 56 kDa subunit contains a region similar to a zinc-binding domain and a region sharing homology with the catalytic loop of a kinase domain. TFIIE binds to RNA polymerase II in solution and joins the preinitiation complex probably concomitant with RNA polymerase II and TFIIF.

α -TFIIE β is an affinity-purified rabbit polyclonal antibody raised against a recombinant full-length p34 subunit of the TFIIE complex.

Specificity

α -TFIIE β reacts with the p34 protein in HeLa nuclear extract (cat.# PR-777) by Western blotting. Recommended dilution range for Western blot analysis: 1:200 - 1:1000

Selected References:

- Ohkuma et al. (1990) *Proc. Natl. Acad. Sci. USA* **87**:9163.
Inostroza et al. (1991) *J. Biol. Chem.* **266**:9304.
Peterson et al. (1991) *Nature* **354**:369.
Flores et al. (1990) *J. Biol. Chem.* **265**:5629.