

α -TFIIA (anti-p55)

anti-p55 subunit of human Transcription Factor II A

rabbit, polyclonal

Cat. No.	Amount
ABD-002	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 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.

α -TFIIA is an affinity-purified rabbit polyclonal antibody raised against a recombinant full-length p55 subunit of the TFIIA complex.

Specificity

α -TFIIA reacts with the p55 protein in HeLa nuclear extract (cat.# PR-777) by Western blotting. Recommended dilution range for Western blot analysis: 1:200 - 1:1000.

Selected References:

- Buratowski et al. (1989) *Cell* **56**:549.
Ranish et al. (1991) *J. Biol. Chem.* **266**:19320.
DeJong et al. (1993) *Genes & Dev.* **7**:2220.
Ma et al. (1993) *Genes & Dev.* **7**:246