

α -RNA Pol II (anti-p33) anti-p33 subunit of RNA Polymerase II rabbit, polyclonal

Cat. No.	Amount
ABD-009	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

hRPB5 (p33) is a highly conserved subunit shared by all three RNA polymerases. It has been shown to be in close contact to promoter DNA when Pol II is recruited into the preinitiation complex. RPB5 has also been implicated in direct protein-protein contacts with transcription factor IIB, RAP30 subunit of transcription factor IIF, and gene-specific modulator proteins, such as the hepatitis B virus transactivator protein X or an inhibitor of Pol II, RMP (RPB5-mediating protein). Therefore, RPB5 is facilitating the communication between the Pol II core and a variety of basal and gene-specific transcription factors. In the Pol II complex, RPB5 interacts with RPB3 and the RPB3-RPB5 interaction is intensified in the presence of three subunits, RPB7, RPB8 and RPB11. RPB5 also makes direct contacts with both RPB1 and RPB2, the two large subunits of the polymerase complex.

α -RNA Pol II (p33) is an affinity-purified rabbit polyclonal antibody raised against a recombinant full-length p33 subunit of the polymerase complex.

Specificity

α -RNA Pol II (p33) reacts with the p33 protein in HeLa nuclear extract (cat.# PR-777) by Western blotting. Recommended dilution range for Western blot analysis: 1:200 - 1:1000.

Selected References:

- Woychik et al. (1990) *Genes Dev.* **4**:313.
- Kim et al. (1997) *Proc. Natl. Acad. Sci. USA* **94**:12268.
- Lin et al. (1997) *J. Biol. Chem.* **272**:7132.
- Wei et al. (2001) *J. Biol. Chem.* **276**:12266.
- Dorjsuren et al. (1998) *Mol. Cell. Biol.* **18**:7546.
- Miyao et al. (1996) *Genes Cells* **1**:843.
- Kimura et al. (2000) *Nucleic Acids Res.* **15**:952.