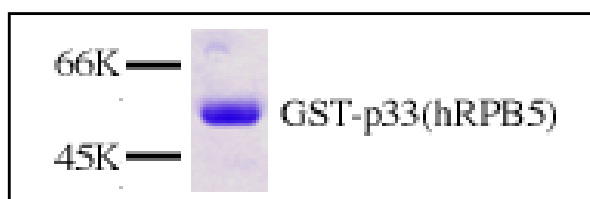


## RNA pol II-hRPB5<sup>GST</sup> RNA Polymerase II, p33 subunit human, recombinant, *E. coli*

Cat. No.	Amount
PR-791	10 µg



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

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 genespecific 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.

Recombinant human RPB5 is isolated from an *E. coli* strain that carries the coding sequence of human RPB5 under the control of a T7 promoter.

RPB5 has been applied in protein-protein interaction assays.

### Selected References:

- Woychik *et al.* (1990) Subunits shared by eukaryotic nuclear RNA polymerases. *Genes Dev.* **4**:313.
- Kim *et al.* (1997) Trajectory of DNA in the RNA polymerase II transcription preinitiation complex. *Proc. Natl. Acad. Sci. USA* **94**:12268.
- Lin *et al.* (1997) Hepatitis B virus X protein is a transcriptional modulator that communicates with transcription factor IIB and the RNA polymerase II subunit 5. *J. Biol. Chem.* **272**:7132.
- Wei *et al.* (2001) Direct interaction between the subunit RAP30 of transcription factor IIF (TFIIF) and RNA polymerase subunit 5, which contributes to the association between TFIIF and RNA polymerase II. *J. Biol. Chem.* **276**:12266.
- Dorjsuren *et al.* (1998) RMP, a novel RNA polymerase II subunit 5-interacting protein, counteracts transactivation by hepatitis B virus X protein. *Mol. Cell. Biol.* **18**:7546.
- Miyao *et al.* (1996) Molecular assembly of RNA polymerase II from the fission yeast *Schizosaccharomyces pombe*: subunit-subunit contact network involving Rpb5. *Genes Cells* **1**:843.
- Kimura *et al.* (2000) Involvement of multiple subunit-subunit contacts in the assembly of RNA polymerase II. *Nucleic Acids Res.* **15**:952.