

## RNA pol II-CTD<sup>GST</sup>

### RNA Polymerase II, Carboxy-terminal Domain

human, recombinant, *E. coli*

Cat. No.	Amount
PR-796	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.

#### Application

GST-CTD has been applied in protein-protein interaction assays.

#### Activity

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

#### Purity

> 95% by SDS-PAGE

#### Description

The carboxy-terminal repeat domain (CTD) of the largest subunit of RNA Pol II contains tandem repeats of a heptapeptide sequence Tyr-Ser-Pro-Thr-Ser-Pro-Ser which is highly conserved among eukaryotic organisms. There are two forms of RNA Pol II *in vivo*, designated IIO, which is extensively phosphorylated at the CTD, and IIA, which is not phosphorylated. The IIA form preferentially enters the pre-initiation complex (PIC), whereas IIO is found in the elongating complex. The kinase activity of TFIIH can mediate CTD phosphorylation, although other kinases, including Cdc2, Ctk1, the Srb10-Srb11 kinase-cyclin pair, and P-TEFβ, have also been implicated in CTD phosphorylation. A phosphatase responsible for the dephosphorylation of the CTD has also been identified. CTD phosphatase activity is regulated by TFIIIB and TFIIIF. The CTD has also been implicated in pre-mRNA processing, most likely functioning as a platform for the recruitment and assembly of factors involved in premRNA processing.

Recombinant GST-CTD is isolated from an *E. coli* strain that carries the coding sequence of human RNA Pol II carboxy-terminal domain under the control of a T7 promoter.

#### Selected References:

- Meinhart *et al.* (2004) Recognition of RNA polymerase II carboxyterminal domain by 3'-RNA-processing factors. *Nature* **430**:223.
- Lu *et al.* (1991) The nonphosphorylated form of RNA polymerase II preferentially associates with the preinitiation complex. *Proc. Natl. Acad. Sci. USA* **88**:10004.
- Dahmus (1994) The role of multisite phosphorylation in the regulation of RNA polymerase II activity. *Prog. Nucleic Acid Res. Mol. Biol.* **48**:143.
- O'Brien *et al.* (1994) Phosphorylation of RNA polymerase II Cterminal domain and transcriptional elongation. *Nature* **370**:75.
- Feaver *et al.* (1991) CTD kinase associated with yeast RNA polymerase II initiation factor b. *Cell* **67**:1223.
- Cisek *et al.* (1989) Phosphorylation of RNA polymerase by the murine homologue of the cell-cycle control protein cdc2. *Nature* **339**:679.
- Lee *et al.* (1991) CTD kinase large subunit is encoded by CTK1, a gene required for normal growth of *Saccharomyces cerevisiae*. *Gene Expr.* **1**:149.
- Liao *et al.* (1995) A kinase-cyclin pair in the RNA polymerase II holoenzyme. *Nature* **374**:193.
- Marshall *et al.* (1996) Control of RNA polymerase II elongation potential by a novel carboxyl-terminal domain kinase. *J. Biol. Chem.*

## RNA pol II-CTD<sup>GST</sup>

### RNA Polymerase II, Carboxy-terminal Domain

#### human, recombinant, *E. coli*

**271:**27176.

Chambers *et al.* (1994) Purification and characterization of a phosphatase from HeLa cells which dephosphorylates the C-terminal domain of RNA polymerase II. *J. Biol. Chem.* **269**:26243.

Chambers *et al.* (1995) The activity of COOH-terminal domain phosphatase is regulated by a docking site on RNA polymerase II and by the general transcription factors IIF and IIB. *J. Biol. Chem.*

**270:**14962.

McCracken *et al.* (1997) The C-terminal domain of RNA polymerase II couples mRNA processing to transcription. *Nature* **385**:357.

Cho *et al.* (2010) Xylocydine, a novel Cdk inhibitor, is an effective inducer of apoptosis in hepatocellular carcinoma cells in vitro and in vivo. *Cancer Letters* **287(2)**:196-206.