

## SV40 T Antigen

### Simian Virus 40 - Large Antigen

human, recombinant, Sf9 insect cells

Cat. No.	Amount
PR-841	5 µg

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

#### Avoid freeze / thaw cycles

#### Form

Liquid. Supplied in 20 mM Tris-HCl pH 8.0, 20% glycerol, 100 mM KCl, 0.2 mM EDTA and 1 mM DTT.

#### Activity

50-500 ng are sufficient for *in vitro* unwinding assay and 100 ng are sufficient for a protein-protein interaction assay.

#### Application

Recombinant SV40 T antigen can be used for

- 1) DNA unwinding assay,
- 2) DNA replication assay,
- 3) DNA binding assays,
- 4) protein-protein interaction assays and
- 5) cell growth assay.

#### Molecular Weight

82 kDa

#### Purity

> 95% by SDS-PAGE

#### Description

The large T antigen (TAg) of Simian virus 40 acts as an initiator of DNA replication, participates in cellular transformation, and induces cell growth. A number of functional domains have been characterized including the N-terminal J domain, nuclear localization signal (NLS), internal helicase domain, Rb and topoisomerase I interacting domain, p53 and ATP binding domain.

The p53-interacting domain is also essential for binding p300/CBP. Phosphorylation of TAg by several kinases may regulate its function both positively and negatively.

#### Selected References:

- Kim *et al.* (2001) Structural basis for the inactivation of retinoblastoma tumor suppressor by SV40 large T antigen. *EMBO J.* **20**:295.
- Chao *et al.* (2000) Loss of p19(ARF) eliminates the requirement for the pRB-binding motif in simian virus 40 large T antigen-mediated transformation. *Mol. Cell. Biol.* **20**:7624.
- Lill *et al.* (1997) Binding and modulation of p53 by p300/CBP coactivators. *Nature* **387**:823.
- Pan *et al.* (1996) Transcription-positive cofactor 4 forms complexes with HSSB (RPA) on single-stranded DNA and influences HSSB-dependent enzymatic synthesis of simian virus 40 DNA. *J. Biol. Chem.* **271**:22111.
- Quartin *et al.* (1994) The amino-terminal functions of the simian virus 40 large T antigen are required to overcome wild-type p53-mediated growth arrest of cells. *J. Virol.* **68**:1334.
- Marton *et al.* (1993) Topoisomerase activity associated with SV40 large tumor antigen. *Nucleic Acids Res.* **21**:1689.
- Moarefi *et al.* (1993) Mutation of the cyclin-dependent kinase phosphorylation site in simian virus 40 (SV40) large T antigen specifically blocks SV40 origin DNA unwinding. *J. Virol.* **67**:4992.
- Fanning (1992) Simian virus 40 large T antigen: the puzzle, the pieces, and the emerging picture. *J. Virol.* **66**:1289.