

α -PC4
anti-Positive Cofactor 4
rabbit, serum

Cat. No.	Amount
ABD-017	100 μ l

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied as serum.

Description

The human PC4 (also called p15) is a non-TAF transcription coactivator that mediates activator-dependent transcription by RNA polymerase II *in vitro* through most tested activators. The function of PC4 is apparently through interactions with transcriptional activators and the basal transcription machinery. It is negatively regulated by casein kinase II phosphorylation both *in vivo* and *in vitro*. PC4 strongly binds single stranded DNA and regulates HSSB (RPA)-dependent SV40 DNA replication. Recent studies indicated that PC4 can be acetylated by several histone acetyltransferase.

α -PC4 is rabbit polyclonal serum raised against the re-combinant full-length PC4 protein.

Specificity

α -PC4 reacts with the PC4 protein in HeLa nuclear extract (cat.# PR-777) by Western blotting. Recommended dilution range for Western blot analysis: 1:2000 - 1:5000.

Selected References:

- Ge et al. (1994) *Cell* **78**:513.
- Ge et al. (1994) *Proc. Natl. Acad. Sci. USA* **91**:12691.
- Kretzschmar et al. (1994) *Cell* **78**:525.
- Wu et al. (1998) *J. Biol. Chem.* **273**:12492.
- Brandsen et al. (1997) *Nat. Struct. Biol.* **4**:900.
- Werten et al. (1998) *EMBO J.* **17**:5103.
- Ge H (2000) *Nucleic Acids Res.* **28**:e3.