

P300/CBP Antibody [#C0290]

- Catalog Number:** C0290-1, C0290-2, C0290
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** Q09472/Q92793
- All Names:** CREB-binding protein, CREBBP
E1A-associated protein p300, EP300, P300 transcription coactivator homolog
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** The antiserum was produced against synthesized peptide derived from human P300/CBP.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** P300/CBP antibody detects endogenous levels of P300/CBP associated factor (P/CAF).
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1: 10000
- References:** Gerd A. Blobel, J. Leukoc. Biol., Apr 2002; 71: 545.
Eric Kalkhoven, Hum. Mol. Genet., Feb 2003; 12: 441 - 450.
Wei-fang Shen, Mol. Cell. Biol., Nov 2001; 21: 7509 - 7522.
Frederick C. Nucifora, Science, Mar 2001; 291: 2423.