

## p300 Antibody [#C0289]

<b>Catalog Number:</b>	C0289
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q09472
<b>Other Names:</b>	E1A-associated protein p300, EP300, P300 transcription coactivator homolog
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from N-terminal of human p300.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	p300 antibody detects endogenous levels of total p300 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
<b>References:</b>	Wuchao Yuan, J. Biol. Chem., Apr 1996; 271: 9009. E. Tory Manning, Mol. Cell. Biol., Jun 2001; 21: 3876 - 3887. Jiwen Li, Mol. Cell. Biol., Mar 2000; 20: 2031 - 2042. Gopalswamy Jayaraman, J. Biol. Chem., Jun 1999; 274: 17342.