

p300 (Ab-89) Antibody [#B0528]

Catalog Number:	B0528
Concentration:	1mg/ml
Swiss-Prot No.:	Q09472
Other Names:	E1A-associated protein p300, EC 2.3.1.48, EP300, P300 transcription coactivator homolog
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg^{2+} and Ca^{2+}), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized non-phosphopeptide derived from human p300 around the phosphorylation site of serine 89 (S-G-S ^P -S-P).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	p300 (Ab-89) antibody detects endogenous levels of total p300 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	P Yaciuk, Mol. Cell. Biol., Apr 1991; 11: 2116 - 2124. JM Routes, J. Immunol., Feb 1996; 156: 1055. Michelle L. Kingsley-Kallesen, J. Biol. Chem., Nov 1999; 274: 34020. Leandro Cerchietti, Blood (ASH Annual Meeting Abstracts), Nov 2006; 108: 1413.