

APEX1 Antibody [#C10493]

Catalog Number:	C10493
Concentration:	1mg/ml
Swiss-Prot No.:	P27695
Other Names:	AP endonuclease 1, APE, APE1, APEN, APEX, APEX nuclease, APEX1, APX, DNA-(apurinic or apyrimidinic site) lyase, EC 4.2.99.18, HAP1, REF-1 protein, REF1
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 peptide derived from internal of human APEX1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	APEX1 antibody detects endogenous levels of total APEX1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Robson C.N., Nucleic Acids Res. 20:4417-4421(1992). Xanthoudakis S., EMBO J. 11:3323-3335(1992). Zhao B., Nucleic Acids Res. 20:4097-4098(1992).