

KAP1 Antibody [#C10527]

Catalog Number:	C10527
Concentration:	1mg/ml
Swiss-Prot No.:	P31321
Other Names:	KAP1, PRKAR1B, cAMP-dependent protein kinase type I-beta regulatory subunit
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 KAP1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	KAP1 antibody detects endogenous levels of total KAP1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Solberg R., Biochem. Biophys. Res. Commun. 176:166-172(1991). Solberg R., Exp. Cell Res. 214:595-605(1994). Zhan X., Anal. Biochem. 354:279-289(2006).