

KAPCB Antibody [#C10453]

Catalog Number:	C10453
Concentration:	1mg/ml
Swiss-Prot No.:	P22694
Other Names:	EC 2.7.11.11, KAPB, KAPCB, PKA C-beta, PKA-beta, cAMP-dependent protein kinase, beta-catalytic 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 KAPCB.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	KAPCB antibody detects endogenous levels of total KAPCB protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Beebe S.J., Mol. Endocrinol. 4:465-475(1990). Gregory S.G., Nature 441:315-321(2006). Wu K.-J., Oncogene 21:7872-7882(2002).