

PRKY Antibody [#C18602]

Catalog Number:	C18602
Concentration:	1mg/ml
Swiss-Prot No.:	O43930
Other Names:	EC 2.7.11.1; Serine/threonine-protein kinase PRKY
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 PRKY.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PRKY antibody detects endogenous levels of total PRKY protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Schiebel K., Hum. Mol. Genet. 6:1985-1989(1997). The MGC Project Team; Genome Res. 14:2121-2127(2004). Schiebel K., Cytogenet. Cell Genet. 76:49-52(1997).