

PRKAG1/2/3 Antibody [#C10728]

Catalog Number:	C10728
Concentration:	1mg/ml
Swiss-Prot No.:	P54619
Other Names:	5'-AMP-activated protein kinase subunit gamma-1, AMPK gamma-1 chain, AMPKg, PRKAG1
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 PRKAG1/2/3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PRKAG1/2/3 antibody detects endogenous levels of total PRKAG1/2/3 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Gao G., J. Biol. Chem. 271:8675-8681(1996). The MGC Project Team, Genome Res. 14:2121-2127(2004). Baba M., Proc. Natl. Acad. Sci. U.S.A. 103:15552-15557(2006).