

PIP5K1C Antibody [#C17692]

Catalog Number:	C17692
Concentration:	1mg/ml
Swiss-Prot No.:	O60331
Other Names:	diphosphoinositide kinase; EC 2.7.1.68; phosphatidylinositol-4-phosphate 5-kinase I-gamma; phosphatidylinositol-4-phosphate 5-kinase, type I, gamma; PI51C; PIP5Kgamma; PIPK I-gamma; PtdIns(4)P-5-kinase; type I PIP kinase
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 PIP5K1C.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PIP5K1C antibody detects endogenous levels of total PIP5K1C protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Nagase T., DNA Res. 5:31-39(1998). Grimwood J., Nature 428:529-535(2004). Di Paolo G., Nature 420:85-89(2002).