

PLCG2(Phospho-Tyr1217) Antibody [#8A0553]

Catalog Number:	8A0553
Concentration:	1mg/ml
Origin of Antibody:	Rabbit
Swiss-Prot No.:	P16885
Other Names:	1-phosphatidylinositol-4,5-bisphosphate phosphodiesterase gamma 2, EC 3.1.4.11, PLC-IV, PLC-gamma2, Phospholipase C-gamma-2
All Sites:	Human: Tyr1217; Mouse: Tyr1217; Rat: Tyr1217
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 phosphopeptide derived from human PLCG2 around the phosphorylation site of tyrosine 1217 (L-F-L-YP-D-T).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.
Specificity:	PLCG2 (Phospho-Tyr1217) antibody detects endogenous levels of PLCG2 only when phosphorylated at tyrosine 1217.
Reactivity:	Human, Mouse, Rat.
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:100~1:500 ELISA: 1:1000
References:	Anouk Caraux, Blood, Feb 2006; 107: 994 - 1002. Yeun Ju Kim, Mol. Cell. Biol., Nov 2004; 24: 9986 - 9999. Jill C. Todt, J. Leukoc. Biol., Apr 2004; 75: 705 - 713. Marie-Pierre Gratacap, J. Biol. Chem., Sep 1998; 273: 24314.