

PLA2G4C Antibody [#C15308]

Catalog Number:	C15308
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UP65
Other Names:	cPLA2-gamma; cytosolic phospholipase A2 gamma precursor; phospholipase A2 group IVC
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 PLA2G4C.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PLA2G4C antibody detects endogenous levels of total PLA2G4C protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Underwood K.W., J. Biol. Chem. 273:21926-21932(1998). Pickard R.T., J. Biol. Chem. 274:8823-8831(1999). Livingston R.J., Submitted (NOV-2003) to the EMBL/GenBank/DDBJ databases.