

AGPAT5 Antibody [#C14006]

Catalog Number:	C14006
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NUQ2
Other Names:	1-acyl-sn-glycerol-3-phosphate acyltransferase epsilon; 1-AGP acyltransferase 5; 1-AGPAT 5; 1-AGPAT5; EC 2.3.1.51; PLCE
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 AGPAT5.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	AGPAT5 antibody detects endogenous levels of total AGPAT5 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Wiemann S., Genome Res. 11:422-435(2001). Ota T., Nat. Genet. 36:40-45(2004). The MGC Project Team; Genome Res. 14:2121-2127(2004).