

PLD3 Antibody [#C17716]

Catalog Number:	C17716
Concentration:	1mg/ml
Swiss-Prot No.:	Q8IV08
Other Names:	Phospholipase D3; EC 3.1.4.4; PLD 3; Choline phosphatase 3; Phosphatidylcholine-hydrolyzing phospholipase D3; HindIII K4L homolog; Hu-K4
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 PLD3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PLD3 antibody detects endogenous levels of total PLD3 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 94%, Positives = 94%); Rat (Identities = 94%, Positives = 94%)
Applications:	IHC: 1:50~1:100 ELISA: 1:5000
References:	Cao J.X., Virus Res. 48:11-18(1997). The MGC Project Team; Genome Res. 14:2121-2127(2004). Munck A., FEBS J. 272:1718-1726(2005).