

LMTK3 Antibody [#C11528]

Catalog Number :	C11528
Concentration:	1mg/ml
Swiss-Prot No.:	Q96Q04
Other Names:	Serine/threonine-protein kinase LMTK3 [Precursor], EC 2.7.11.1, Lemur tyrosine kinase 3, LMTK3, KIAA1883, TYKLM3
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 LMTK3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	LMTK3 antibody detects endogenous levels of total LMTK3 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Nagase T., DNA Res. 8:179-187(2001). The MGC Project Team, Genome Res. 14:2121-2127(2004).