

LDOCL Antibody [#C11219]

Catalog Number:	C11219
Concentration:	1mg/ml
Swiss-Prot No.:	Q6ICC9
Other Names:	Protein LDOC1L, Leucine zipper protein down-regulated in cancer cells-like, Mammalian retrotransposon-derived protein 6, LDOC1L, MAR6, MART6
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 LDOCL.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	LDOCL antibody detects endogenous levels of total LDOCL protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Collins J.E., Genome Biol. 5:RESEARCH84.1-RESEARCH84.11(2004). Dunham I., Nature 402:489-495(1999). Ota T., Nat. Genet. 36:40-45(2004).