

MELK Antibody [#C11043]

Catalog Number:	C11043
Concentration:	1mg/ml
Swiss-Prot No.:	Q14680
Other Names:	Maternal embryonic leucine zipper kinase, EC 2.7.11.1, hMELK, Protein kinase PK38, hPK38, KIAA0175
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 MELK.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MELK antibody detects endogenous levels of total MELK protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Nagase T., DNA Res. 3:17-24(1996). Seong H.-A., Biochem. J. 361:597-604(2002). Vulsteke V., J. Biol. Chem. 279:8642-8647(2004).