

MAEA Antibody [#C11253]

Catalog Number:	C11253
Concentration:	1mg/ml
Swiss-Prot No.:	Q7L5Y9
Other Names:	Macrophage erythroblast attacher, Erythroblast macrophage protein, Cell proliferation-inducing gene 5 protein, Lung cancer-related protein 10, HLC-10, MAEA, EMP, HLC10, PIG5
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 MAEA.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MAEA antibody detects endogenous levels of total MAEA protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Ota T., Nat. Genet. 36:40-45(2004). Hanspal M., Blood 92:2940-2950(1998). Bala S., Biochem. Biophys. Res. Commun. 342:1040-1048(2006).