

NOL10 Antibody [#C17154]

Catalog Number:	C17154
Concentration:	1mg/ml
Swiss-Prot No.:	Q9BSC4
Other Names:	FLJ13938; FLJ14075
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 C-terminal of human NOL10.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	NOL10 antibody detects endogenous levels of total NOL10 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Ota T., Nat. Genet. 36:40-45(2004). Hillier L.W., Nature 434:724-731(2005). The MGC Project Team; Genome Res. 14:2121-2127(2004).