

CALML5 Antibody [#C14863]

Catalog Number:	C14863
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NZT1
Other Names:	Calmodulin-like protein 5; Calmodulin-like skin protein
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 CALML5.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CALML5 antibody detects endogenous levels of total CALML5 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:20000
References:	Mehul B., J. Biol. Chem. 275:12841-12847(2000). The MGC Project Team; Genome Res. 14:2121-2127(2004).