

## PKCB1 Antibody [#C11882]

<b>Catalog Number:</b>	C11882
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9ULU4
<b>Other Names:</b>	CTCL tumor antigen se14-3, Cutaneous T-cell lymphoma associated antigen se14-3, KIAA1125, PKCB, PKCB1, PRKCBP1, Protein kinase C binding protein 1, ZMYND8, Zinc finger MYND domain containing protein 8
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human PKCB1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	PKCB1 antibody detects endogenous levels of total PKCB1 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Fossey S.C., Mamm. Genome 11:919-925(2000). Eichmueller S., Proc. Natl. Acad. Sci. U.S.A. 98:629-634(2001). Deloukas P., Nature 414:865-871(2001).