

## PDLIM1 Antibody [#C10012]

<b>Catalog Number:</b>	C10012
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	O00151
<b>Other Names:</b>	C-terminal LIM domain protein 1, CLP36, Elfin, LIM domain protein CLP-36, PDL1, PDLI1, PDLIM1, PDZ and LIM domain protein 1
<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 PDLIM1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	PDLIM1 antibody detects endogenous levels of total PDLIM1 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Kotaka M., J. Cell. Biochem. 72:279-285(1999). Bauer K., Blood 96:4236-4245(2000). Deloukas P., Nature 429:375-381(2004).