

## CBLN1 Antibody [#C15056]

<b>Catalog Number:</b>	C15056
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P23435
<b>Other Names:</b>	Cerebellin-1; Precerebellin
<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 CBLN1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	CBLN1 antibody detects endogenous levels of total CBLN1 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000   ELISA: 1:40000
<b>References:</b>	Urade Y., Proc. Natl. Acad. Sci. U.S.A. 88:1069-1073(1991). Ota T., Nat. Genet. 36:40-45(2004). Mural R.J., Submitted (JUL-2005) to the EMBL/GenBank/DDBJ databases.