

APLP2 Antibody [#C14458]

Catalog Number:	C14458
Concentration:	1mg/ml
Swiss-Prot No.:	Q06481
Other Names:	Amyloid protein homolog; Amyloid-like protein 2 precursor; APLP2; APP2; APPH; APPL2; CDEBP; CDEI-box binding 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 APLP2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	APLP2 antibody detects endogenous levels of total APLP2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%) Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Sprecher C.A., Biochemistry 32:4481-4486(1993). Wasco W., Nat. Genet. 5:95-99(1993). von der Kammer H., DNA Cell Biol. 13:1137-1143(1994).