

CD2L1 Antibody [#C10442]

Catalog Number:	C10442
Concentration:	1mg/ml
Swiss-Prot No.:	P21127
Other Names:	CD2L1, CDC2L1, CDK11, CLK-1, Cell division cycle 2-like 1, EC 2.7.11.22, Galactosyltransferase associated protein kinase p58/GTA, PITSLREA, PK58, p58 CLK-1
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 N-terminal of human CD2L1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CD2L1 antibody detects endogenous levels of total CD2L1 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Bunnell B.A., Proc. Natl. Acad. Sci. U.S.A. 87:7467-7471(1990). Eipers P.G., Genomics 13:613-621(1992). Xiang J., J. Biol. Chem. 269:15786-15794(1994).