

60S Ribosomal Protein L10 Antibody [#C0391]

Catalog Number:	C0391
Concentration:	1mg/ml
Swiss-Prot No.:	P27635
Other Names:	60S ribosomal protein L10; Protein QM; Tumor suppressor QM; Laminin receptor homolog; RPL10; DXS648E, QM
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 60S Ribosomal Protein L10.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	60S Ribosomal Protein L10 antibody detects endogenous levels of total 60S Ribosomal Protein L10 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Dowdy S.F., Nucleic Acids Res. 19:5763-5769(1991). van den Ouweland A.M.W., Hum. Mol. Genet. 1:269-273(1992). Kaneko K., Hum. Mol. Genet. 1:529-533(1992).