

RPL18 Antibody [#C14154]

Catalog Number:	C14154
Concentration:	1mg/ml
Swiss-Prot No.:	Q07020
Other Names:	ribosomal protein L18; RL18
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 RPL18.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RPL18 antibody detects endogenous levels of total RPL18 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:	Puder M., Biochim. Biophys. Acta 1216:134-136(1993). Yoshihama M., Genome Res. 12:379-390(2002). Odintsova T.I., J. Protein Chem. 22:249-258(2003).