

RS27L Antibody [#C14115]

Catalog Number:	C14115
Concentration:	1mg/ml
Swiss-Prot No.:	Q71UM5
Other Names:	40S ribosomal protein S27-like protein; RPS27L
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 RS27L.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RS27L antibody detects endogenous levels of total RS27L protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Zhang Q.-H., Genome Res. 10:1546-1560(2000). The MGC Project Team; Genome Res. 14:2121-2127(2004).