

RPS11 Antibody [#C14094]

Catalog Number:	C14094
Concentration:	1mg/ml
Swiss-Prot No.:	P62280
Other Names:	40S ribosomal protein S11; ribosomal protein S11; RS11
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 RPS11.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RPS11 antibody detects endogenous levels of total RPS11 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Lott J.B., Nucleic Acids Res. 16:1205-1205(1988). Higa S., Gene 240:371-377(1999). The MGC Project Team; Genome Res. 14:2121-2127(2004).