

RPS8 Antibody [#C14126]

Catalog Number:	C14126
Concentration:	1mg/ml
Swiss-Prot No.:	P62241
Other Names:	40S ribosomal protein S8; OK/SW-cl.83; ribosomal protein S8; RS8
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 RPS8.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RPS8 antibody detects endogenous levels of total RPS8 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Davies B., Genomics 15:68-75(1993). Shichijo S., Submitted (MAY-2001) to the EMBL/GenBank/DDBJ databases. The MGC Project Team; Genome Res. 14:2121-2127(2004).