

RPS13 Antibody [#C14096]

Catalog Number:	C14096
Concentration:	1mg/ml
Swiss-Prot No.:	P62277
Other Names:	40S ribosomal protein S13
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 RPS13.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RPS13 antibody detects endogenous levels of total RPS13 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Rush J., Nat. Biotechnol. 23:94-101(2005). Olsen J.V., Cell 127:635-648(2006). Kenmochi N., Biochem. Biophys. Res. Commun. 228:371-374(1996).