

RPS3 Antibody [#C14118]

Catalog Number:	C14118
Concentration:	1mg/ml
Swiss-Prot No.:	P23396
Other Names:	40S ribosomal protein S3; RS3
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 RPS3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RPS3 antibody detects endogenous levels of total RPS3 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
References:	Yoshihama M., Genome Res. 12:379-390(2002). Nousiainen M., Proc. Natl. Acad. Sci. U.S.A. 103:5391-5396(2006). Molina H., Proc. Natl. Acad. Sci. U.S.A. 104:2199-2204(2007).