

RPS21 Antibody [#C14107]

Catalog Number:	C14107
Concentration:	1mg/ml
Swiss-Prot No.:	P63220
Other Names:	40S ribosomal protein S21
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 RPS21.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RPS21 antibody detects endogenous levels of total RPS21 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:	Yoshihama M., Genome Res. 12:379-390(2002). Deloukas P., Nature 414:865-871(2001). Kenmochi N., Genome Res. 8:509-523(1998).