

MRPL47 Antibody [#C14083]

Catalog Number:	C14083
Concentration:	1mg/ml
Swiss-Prot No.:	Q9HD33
Other Names:	39S ribosomal protein L47, mitochondrial; nasopharyngeal carcinoma metastasis-related 1; NCM1; RM47
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 MRPL47.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MRPL47 antibody detects endogenous levels of total MRPL47 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Muzny D.M., Nature 440:1194-1198(2006). Mural R.J., Submitted (SEP-2005) to the EMBL/GenBank/DDBJ databases. The MGC Project Team; Genome Res. 14:2121-2127(2004).