

## MRPL12 Antibody [#C14056]

<b>Catalog Number:</b>	C14056
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P52815
<b>Other Names:</b>	39S ribosomal protein L12, mitochondrial; 5c5-2; L12mt; MRP-L12; RM12; RPML12
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human MRPL12.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	MRPL12 antibody detects endogenous levels of total MRPL12 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%) Mouse (Identities = 100%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Marty L., J. Biol. Chem. 271:11468-11476(1996). Liu J., Submitted (APR-1998) to the EMBL/GenBank/DDBJ databases. The MGC Project Team; Genome Res. 14:2121-2127(2004).