

MRPL22 Antibody [#C14067]

Catalog Number:	C14067
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NWU5
Other Names:	39S ribosomal protein L22, mitochondrial; L22mt; mitochondrial ribosomal protein L22; MRP-L22; MRP-L25; RM22; RPML25
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 MRPL22.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MRPL22 antibody detects endogenous levels of total MRPL22 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Zhang Q.-H., Genome Res. 10:1546-1560(2000). Ota T., Nat. Genet. 36:40-45(2004). The German cDNA consortium; Submitted (JUN-2005) to the EMBL/GenBank/DDBJ databases.