

MRPL46 Antibody [#C14082]

Catalog Number:	C14082
Concentration:	1mg/ml
Swiss-Prot No.:	Q9H2W6
Other Names:	mitochondrial ribosomal protein l46
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 MRPL46.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MRPL46 antibody detects endogenous levels of total MRPL46 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 82%, Positives = 94%); Rat (Identities = 82%, Positives = 94%)
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
References:	Carim-Todd L., DNA Seq. 12:91-96(2001). The MGC Project Team; Genome Res. 14:2121-2127(2004). Koc E.C., J. Biol. Chem. 276:43958-43969(2001).