

MRPL21 Antibody [#C14066]

Catalog Number:	C14066
Concentration:	1mg/ml
Swiss-Prot No.:	Q7Z2W9
Other Names:	39S ribosomal protein L21, mitochondrial; L21mt; RP-L21
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 MRPL21.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MRPL21 antibody detects endogenous levels of total MRPL21 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 82%, Positives = 94%)
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
References:	The MGC Project Team; Genome Res. 14:2121-2127(2004). Colinge J., Submitted (OCT-2008) to UniProtKB.