

MRPL11 Antibody [#C14055]

Catalog Number:	C14055
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y3B7
Other Names:	cgi-113; cgi-113 protein; l11mt; mgc111024; mitochondrial ribosomal protein l11; mitochondrial ribosomal protein l11 (l11mt); mitochondrial ribosomal protein l11 isoform a; mitochondrial ribosomal protein l11 isoform a variant; mitochondrial ribosomal protein l11 isoform b; mrp-l11; mrpl11; rm11
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 MRPL11.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MRPL11 antibody detects endogenous levels of total MRPL11 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
References:	Suzuki T., J. Biol. Chem. 276:21724-21736(2001). Lai C.-H., Genome Res. 10:703-713(2000). The MGC Project Team; Genome Res. 14:2121-2127(2004).