

MRPL50 Antibody [#C16675]

Catalog Number:	C16675
Concentration:	1mg/ml
Swiss-Prot No.:	Q8N5N7
Other Names:	flj20493; flj21990; mitochondrial ribosomal protein l50; mrp-l50; mrpl50; rm50
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 MRPL50.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MRPL50 antibody detects endogenous levels of total MRPL50 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Ota T., Nat. Genet. 36:40-45(2004). Humphray S.J., Nature 429:369-374(2004). The MGC Project Team; Genome Res. 14:2121-2127(2004).