

MRPL44 Antibody [#C14080]

Catalog Number:	C14080
Concentration:	1mg/ml
Swiss-Prot No.:	Q9H9J2
Other Names:	39S ribosomal protein L44, mitochondrial; EC 3.1.26.-; RM44
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 MRPL44.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MRPL44 antibody detects endogenous levels of total MRPL44 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
References:	Ota T., Nat. Genet. 36:40-45(2004). Ebert L., Submitted (JUN-2004) to the EMBL/GenBank/DDBJ databases. Hillier L.W., Nature 434:724-731(2005).