

MRPS25 Antibody [#C16655]

Catalog Number:	C16655
Concentration:	1mg/ml
Swiss-Prot No.:	P82663
Other Names:	mitochondrial 28S ribosomal protein S25; mitochondrial ribosomal protein S25; MRP-S25; RPMS25; RT25; S25mt
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 MRPS25.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MRPS25 antibody detects endogenous levels of total MRPS25 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 92%); Rat (Identities = 92%, Positives = 92%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Ota T., Nat. Genet. 36-40-45(2004). Hattori A., DNA Res. 7:357-366(2000). Koc E.C., J. Biol. Chem. 275:32585-32591(2000).