

MRPS16 Antibody [#C14035]

Catalog Number:	C14035
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y3D3
Other Names:	mitochondrial ribosomal protein S16; MRP-S16; RPMS16; RT16; S16mt
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 MRPS16.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MRPS16 antibody detects endogenous levels of total MRPS16 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
References:	Lai C.-H., Genome Res. 10:703-713(2000). Suzuki T., J. Biol. Chem. 276:33181-33195(2001). Kenmochi N., Genomics 77:65-70(2001).