

MRPS9 Antibody [#C14045]

Catalog Number:	C14045
Concentration:	1mg/ml
Swiss-Prot No.:	P82933
Other Names:	28S ribosomal protein S9, mitochondrial; MRP-S9; RPMS9; RT09; S9mt
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 MRPS9.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MRPS9 antibody detects endogenous levels of total MRPS9 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%) Mouse (Identities = 98%, Positives = 194%); Rat (Identities = 94%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	The MGC Project Team; Genome Res. 14:2121-2127(2004). Koc E.C., J. Biol. Chem. 276:19363-19374(2001).