

MRPS18C Antibody [#C14039]

Catalog Number:	C14039
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y3D5
Other Names:	28S ribosomal protein S18c, mitochondrial; MRP-S18-1; MRP-S18-c; RT18C
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 MRPS18C.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MRPS18C antibody detects endogenous levels of total MRPS18C protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
References:	Lai C.-H., Genome Res. 10:703-713(2000). The MGC Project Team; Genome Res. 14:2121-2127(2004). Koc E.C., J. Biol. Chem. 276:19363-19374(2001).