

## MRPL20 Antibody [#C14065]

<b>Catalog Number:</b>	C14065
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9BYC9
<b>Other Names:</b>	39S ribosomal protein L20, mitochondrial [Precursor]; L20mt; MRP-L20; MRPL20
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from C-terminal of human MRPL20.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	MRPL20 antibody detects endogenous levels of total MRPL20 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Suzuki T., J. Biol. Chem. 276:21724-21736(2001). Gregory S.G., Nature 441:315-321(2006). The MGC Project Team; Genome Res. 14:2121-2127(2004).