

## RPL10L Antibody [#C14146]

<b>Catalog Number:</b>	C14146
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q96L21
<b>Other Names:</b>	60S ribosomal protein L10-like; RL10L
<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 internal of human RPL10L.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	RPL10L antibody detects endogenous levels of total RPL10L protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Uechi T., Nucleic Acids Res. 30:5369-5375(2002). The MGC Project Team; Genome Res. 14:2121-2127(2004).