

## RPL39 Antibody [#C14181]

<b>Catalog Number:</b>	C14181
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P62891
<b>Other Names:</b>	60S ribosomal protein L39; RPL39
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), 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 RPL39.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	RPL39 antibody detects endogenous levels of total RPL39 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:60000
<b>References:</b>	Tsui S.K.W., Biochem. Mol. Biol. Int. 40:611-616(1996). Otsuka S., Biochim. Biophys. Acta 1308:119-121(1996). Yoshihama M., Genome Res. 12:379-390(2002).