

## RPL28 Antibody [#C14166]

<b>Catalog Number:</b>	C14166
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P46779
<b>Other Names:</b>	FLJ43307; ribosomal protein L28; RL28
<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 RPL28.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	RPL28 antibody detects endogenous levels of total RPL28 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Frigerio J.-M., Biochim. Biophys. Acta 1262:64-68(1995). Ebert L., Submitted (JUN-2004) to the EMBL/GenBank/DDBJ databases. The MGC Project Team; Genome Res. 14:2121-2127(2004).