

## RPL12 Antibody [#C14148]

|                           |                                                                                                                                                   |
|---------------------------|---------------------------------------------------------------------------------------------------------------------------------------------------|
| <b>Catalog Number:</b>    | C14148                                                                                                                                            |
| <b>Concentration:</b>     | 1mg/ml                                                                                                                                            |
| <b>Swiss-Prot No.:</b>    | P30050                                                                                                                                            |
| <b>Other Names:</b>       | RL12; RPL12                                                                                                                                       |
| <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 RPL12.                                                      |
| <b>Purification:</b>      | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.                             |
| <b>Specificity:</b>       | RPL12 antibody detects endogenous levels of total RPL12 protein.                                                                                  |
| <b>Reactivity:</b>        | Human, Mouse                                                                                                                                      |
| <b>Applications:</b>      | WB: 1:500~1:1000   ELISA: 1:40000                                                                                                                 |
| <b>References:</b>        | Chu W., Nucleic Acids Res. 21:749-749(1993).<br>Humphray S.J., Nature 429:369-374(2004).<br>The MGC Project Team; Genome Res. 14:2121-2127(2004). |