

RPL15 Antibody [#C14152]

| | |
|---------------------------|--|
| Catalog Number: | C14152 |
| Concentration: | 1mg/ml |
| Swiss-Prot No.: | P61313 |
| Other Names: | 60S ribosomal protein L15; EC45; RL15 |
| Storage/Stability: | Store at -20°C/1 year |
| Form of Antibody: | Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol. |
| Immunogen: | The antiserum was produced against synthesized peptide derived from internal of human RPL15. |
| Purification: | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen. |
| Specificity: | RPL15 antibody detects endogenous levels of total RPL15 protein. |
| Reactivity: | Human, Mouse, Rat |
| Applications: | WB: 1:500~1:1000 ELISA: 1:40000 |
| References: | Herzog H.; Submitted (NOV-1993) to the EMBL/GenBank/DDBJ databases. Wang Q., Submitted (JUN-2000) to the EMBL/GenBank/DDBJ databases. Yoshihama M., Submitted (MAY-2001) to the EMBL/GenBank/DDBJ databases. |