

RRS1 Antibody [#C18435]

| | |
|---------------------------|--|
| Catalog Number: | C18435 |
| Concentration: | 1mg/ml |
| Swiss-Prot No.: | Q15050 |
| Other Names: | ribosome biogenesis regulatory protein; RRR |
| 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 RRS1. |
| Purification: | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen. |
| Specificity: | RRS1 antibody detects endogenous levels of total RRS1 protein. |
| Reactivity: | Human, Mouse, Rat |
| Applications: | WB: 1:500~1:1000 ELISA: 1:10000 |
| References: | Nagase T., DNA Res. 2:37-43(1995). The MGC Project Team; Genome Res. 14:2121-2127(2004). Bienvenut W.V.; Submitted (AUG-2005) to UniProtKB. |