

RAB7L1 Antibody [#C18271]

| | |
|---------------------------|--|
| Catalog Number: | C18271 |
| Concentration: | 1mg/ml |
| Swiss-Prot No.: | O14966 |
| Other Names: | RAB7, member RAS oncogene family-like 1; RAB7L |
| 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 RAB7L1. |
| Purification: | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen. |
| Specificity: | RAB7L1 antibody detects endogenous levels of total RAB7L1 protein. |
| Reactivity: | Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%) |
| Applications: | WB: 1:500~1:1000 ELISA: 1:20000 |
| References: | Shimizu F., Cytogenet. Cell Genet. 77:261-263(1997). The MGC Project Team, Genome Res. 14:2121-2127(2004). |