

KAP2 Antibody [#C10363]

| | |
|---------------------------|--|
| Catalog Number: | C10363 |
| Concentration: | 1mg/ml |
| Swiss-Prot No.: | P13861 |
| Other Names: | KAP2, MGC3606, PKR2, PRKAR2 |
| 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 KAP2. |
| Purification: | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen. |
| Specificity: | KAP2 antibody detects endogenous levels of total KAP2 protein. |
| Reactivity: | Human |
| Applications: | WB: 1:500~1:1000 ELISA: 1:10000 |
| References: | Oyen O., FEBS Lett. 246:57-64(1989). Foss K.B., Biochim. Biophys. Acta 1350:98-108(1997). Schillace R.V., J. Immunol. 168:1590-1599(2002). |