

TRIP4 Antibody [#C11094]

| | |
|---------------------------|--|
| Catalog Number: | C11094 |
| Concentration: | 1mg/ml |
| Swiss-Prot No.: | Q15650 |
| Other Names: | Activating signal cointegrator 1, ASC-1, Thyroid receptor-interacting protein 4, TRIP-4, TRIP4 |
| 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 TRIP4. |
| Purification: | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen. |
| Specificity: | TRIP4 antibody detects endogenous levels of total TRIP4 protein. |
| Reactivity: | Human, Mouse |
| Applications: | WB: 1:500~1:1000 ELISA: 1:5000 |
| References: | Kim H.J., Mol. Cell. Biol. 19:6323-6332(1999). Lee J.W., Mol. Endocrinol. 9:243-254(1995). Jung D.-J., Mol. Cell. Biol. 22:5203-5211(2002). |