

TRIM24 Antibody [#C10073]

| | |
|---------------------------|--|
| Catalog Number: | C10073 |
| Concentration: | 1mg/ml |
| Swiss-Prot No.: | O15164 |
| Other Names: | TIF1, TIF1A, TRIM24, Transcription intermediary factor 1-alpha, Tripartite motif protein 24, kinase TIF1-alpha |
| 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 C-terminal of human TRIM24. |
| Purification: | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen. |
| Specificity: | TRIM24 antibody detects endogenous levels of total TRIM24 protein. |
| Reactivity: | Human, Mouse |
| Applications: | WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000 |
| References: | Thenot S., J. Biol. Chem. 272:12062-12068(1997). Venturini L., Oncogene 18:1209-1217(1999). Hellal-Levy C., Mol. Endocrinol. 14:1210-1221(2000). |