

ZNF174 Antibody [#C11098]

| | |
|---------------------------|---|
| Catalog Number: | C11098 |
| Concentration: | 1mg/ml |
| Swiss-Prot No.: | Q15697 |
| Other Names: | Zinc finger protein 174, AW-1, Zinc finger and SCAN domain-containing protein 8, ZNF174, ZSCAN8 |
| 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 ZNF174. |
| Purification: | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen. |
| Specificity: | ZNF174 antibody detects endogenous levels of total ZNF174 protein. |
| Reactivity: | Human |
| Applications: | WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000 |
| References: | Williams A.J., J. Biol. Chem. 270:22143-22152(1995). The MGC Project Team, Genome Res. 14:2121-2127(2004). Rush J., Nat. Biotechnol. 23:94-101(2005). |