

TNFA Antibody [#C10265]

| | |
|---------------------------|--|
| Catalog Number: | C10265 |
| Concentration: | 1mg/ml |
| Swiss-Prot No.: | P01375 |
| Other Names: | Protein name Tumor necrosis factor [Precursor], TNF-alpha, Tumor necrosis factor ligand superfamily member 2, TNF-a, Cachectin, Tumor necrosis factor, membrane form, Tumor necrosis factor, soluble form, TNF, TNFA, TNFSF2 |
| 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 TNFA. |
| Purification: | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen. |
| Specificity: | TNFA antibody detects endogenous levels of total TNFA protein. |
| Reactivity: | Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%) |
| Applications: | IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000 |
| References: | Neville M.J., J. Immunol. 162:4745-4754(1999). Xie T., Genome Res. 13:2621-2636(2003). Kim Y.J., Hum. Mol. Genet. 12:2541-2546(2003). |