

FAS Antibody [#C0355]

| | |
|---------------------------|---|
| Catalog Number: | C0355 |
| Concentration: | 1mg/ml |
| Swiss-Prot No.: | P25445 |
| Other Names: | Apo-1, CD95, FasL receptor, TNFRSF6, TNFR6, apoptosis-mediating surface antigen FAS, tumor necrosis factor receptor superfamily member 6 |
| 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 Fas. |
| Purification: | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen. |
| Specificity: | Fas antibody detects endogenous levels of total Fas protein. |
| Reactivity: | Human |
| Applications: | WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000 |
| References: | Michael J. Pinkoski, Am J Physiol Gastrointest Liver Physiol, Mar 2000; 278: 354. Morgan Jenkins, J. Immunol., Aug 1999; 163: 1195. Nicholas Mitsiades, J Natl Cancer Inst, Oct 1999; 91: 1678 - 1684. JOSÉ E. CARDIER, FASEB J, Nov 1999; 13: 1950. |