

hnRNP A2/B1 Antibody [#C10452]

| | |
|---------------------------|------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Catalog Number: | C10452 |
| Concentration: | 1mg/ml |
| Swiss-Prot No.: | P22626 |
| Other Names: | HNRPA2B1, Heterogeneous nuclear ribonucleoproteins A2/B1, ROA2, hnRNP A2 / hnRNP B1 |
| 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 N-terminal of human hnRNP A2/B1. |
| Purification: | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen. |
| Specificity: | hnRNP A2/B1 antibody detects endogenous levels of total hnRNP A2/B1 protein. |
| Reactivity: | Human, Mouse |
| Applications: | WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000 |
| References: | Burd C.G., Proc. Natl. Acad. Sci. U.S.A. 86:9788-9792(1989). Kozu T., Genomics 25:365-371(1995). Biamonti G., Nucleic Acids Res. 22:1996-2002(1994). |