

NFYA Antibody [#C10461]

Catalog Number:	C10461
Concentration:	1mg/ml
Swiss-Prot No.:	P23511
Other Names:	CAAT-box DNA binding protein subunit A, CBF-A, CBF-B, CCAAT-binding transcription factor subunit B, HAP2, HAP2 CCAAT-binding protein, NF-Y protein chain A, NFYA, nuclear transcription factor Y, alpha, transcription factor NF-Y, A subunit
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 NFYA.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	NFYA antibody detects endogenous levels of total NFYA protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Li X.-Y., Nucleic Acids Res. 20:1087-1091(1992). Mungall A.J., Nature 425:805-811(2003). Becker D.M., Proc. Natl. Acad. Sci. U.S.A. 88:1968-1972(1991).