

ELF1 Antibody [#C10538]

Catalog Number:	C10538
Concentration:	1mg/ml
Swiss-Prot No.:	P32519
Other Names:	E74-like factor 1, ELF1, ETS-related transcription factor Elf-1
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 ELF1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ELF1 antibody detects endogenous levels of total ELF1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
References:	Leiden J.M., J. Virol. 66:5890-5897(1992). Thompson C.B., Mol. Cell. Biol. 12:1043-1053(1992). Wang C.Y., Science 260:1330-1335(1993).