

CSTF1 Antibody [#C10882]

Catalog Number:	C10882
Concentration:	1mg/ml
Swiss-Prot No.:	Q05048
Other Names:	CF-1 50 kDa subunit, CSTF 50 kDa subunit, CSTF1, Cleavage stimulation factor, 50 kDa 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 N-terminal of human CSTF1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CSTF1 antibody detects endogenous levels of total CSTF1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Takagaki Y., Proc. Natl. Acad. Sci. U.S.A. 89:1403-1407(1992). Deloukas P., Nature 414:865-871(2001). Kleiman F.E., Science 285:1576-1579(1999).