

CELF-1 Antibody [#C11440]

Catalog Number:	C11440
Concentration:	1mg/ml
Swiss-Prot No.:	Q92879
Other Names:	CELF-1, Bruno-like protein 2, RNA-binding protein BRUNOL-2, CUG triplet repeat RNA-binding protein 1, CUG-BP1, Deadenylation factor CUG-BP, 50 kDa nuclear polyadenylated RNA-binding protein, Embryo deadenylation element-binding protein homolog, EDEN-BP homolog
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 CELF-1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CELF-1 antibody detects endogenous levels of total CELF-1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Timchenko L.T., Nucleic Acids Res. 24:4407-4414(1996). Good P.J., J. Biol. Chem. 275:28583-28592(2000). Takahashi N., J. Biochem. 130:581-587(2001).