

FUBP3 Antibody [#C15743]

Catalog Number:	C15743
Concentration:	1mg/ml
Swiss-Prot No.:	Q96I24
Other Names:	far upstream element-binding protein 3; FUBP3; FUSE binding protein 3
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 FUBP3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	FUBP3 antibody detects endogenous levels of total FUBP3 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Davis-Smyth T., J. Biol. Chem. 271:31679-31687(1996). The MGC Project Team; Genome Res. 14:2121-2127(2004). Bienvenut W.V., Submitted (MAR-2008) to UniProtKB.