

TAF1 Antibody [#C10445]

Catalog Number:	C10445
Concentration:	1mg/ml
Swiss-Prot No.:	P21675
Other Names:	BA2R, CCG1, Cell cycle gene 1 protein, EC 2.7.11.1, P250, T2D1, TAF2A, TAFII-250, TAFII250, TBP-associated factor 250 kDa, Transcription initiation factor TFIID 250 kDa subunit, kinase TAF1
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 TAF1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TAF1 antibody detects endogenous levels of total TAF1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:20000
References:	Sekiguchi T., Mol. Cell. Biol. 11:3317-3325(1991). Ruppert S., Nature 362:175-179(1993). Nolte D., Proc. Natl. Acad. Sci. U.S.A. 100:10347-10352(2003).