

TF3C2 Antibody [#C11403]

Catalog Number:	C11403
Concentration:	1mg/ml
Swiss-Prot No.:	Q8WUA4
Other Names:	General transcription factor 3C polypeptide 2, Transcription factor III C subunit beta, TF3C-beta, TFIIC 110 kDa subunit, TFIIC110, GTF3C2, KIAA0011
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 TF3C2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TF3C2 antibody detects endogenous levels of total TF3C2 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Sinn E., Genes Dev. 9:675-685(1995). Nomura N., DNA Res. 1:27-35(1994). Olsen J.V., Cell 127:635-648(2006).