

TF3C3 Antibody [#C11971]

Catalog Number:	C11971
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y5Q9
Other Names:	TF3C-gamma, TF3C3, TFIIC 102 kDa subunit, TFIIC102, TFIICgamma, TFIIIC2-102, eneral transcription factor 3C polypeptide 3, general transcription factor IIC, polypeptide 3 (102kD), general transcription factor IIC, polypeptide 3, 102kDa, transcription factor IIC, 102 kD, transcription factor IIC-gamma 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 internal of human TF3C3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TF3C3 antibody detects endogenous levels of total TF3C3 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Ota T., Nat. Genet. 36:40-45(2004). The German cDNA consortium; Submitted (JAN-2000) to the EMBL/GenBank/DDBJ databases. Hsieh Y.-J., Mol. Cell. Biol. 19:4944-4952(1999).