

TCF3 Antibody [#C10385]

Catalog Number:	C10385
Concentration:	1mg/ml
Swiss-Prot No.:	P15923
Other Names:	E12, ITF1, Immunoglobulin enhancer binding factor E12/E47, Immunoglobulin transcription factor-1, Kappa-E2-binding factor, TCF-3, TCF3, TFE2, Transcription factor E2-alpha, Transcription factor ITF-1, Transcription factor-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 C-terminal of human TCF3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TCF3 antibody detects endogenous levels of total TCF3 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
References:	Kamps M.P., Cell 60:547-555(1990). Nourse J., Cell 60:535-545(1990). Grimwood J., Nature 428:529-535(2004).