

TCFL5 Antibody [#C11878]

Catalog Number:	C11878
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UL49
Other Names:	Transcription factor-like 5 protein, Cha transcription factor, HPV-16 E2-binding protein 1, E2BP-1, TCFL5, CHA, E2BP1
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 TCFL5.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TCFL5 antibody detects endogenous levels of total TCFL5 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Girones N., J. Clin. Invest. 107:985-993(2001). Deloukas P., Nature 414:865-871(2001). Maruyama O., Cytogenet. Cell Genet. 82:41-45(1998).