

## TLE4 Antibody [#C10876]

<b>Catalog Number:</b>	C10876
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q04727
<b>Other Names:</b>	TIM, TIM1, timeless (Drosophila) homolog
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human TLE4.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	TLE4 antibody detects endogenous levels of total TLE4 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000   ELISA: 1:20000
<b>References:</b>	Nagase T., DNA Res. 6:337-345(1999). Humphray S.J., Nature 429:369-374(2004). Stifani S., Nat. Genet. 2:119-127(1992).