

## RBM26 Antibody [#C11188]

<b>Catalog Number:</b>	C11188
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q5T8P6
<b>Other Names:</b>	RNA-binding motif protein 26, RNA-binding protein 26, CTCL tumor antigen se70-2, C13orf10, PRO1777
<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 C-terminal of human RBM26.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	RBM26 antibody detects endogenous levels of total RBM26 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000   ELISA: 1:10000
<b>References:</b>	Ota T., Nat. Genet. 36:40-45(2004). Dunham A., Nature 428:522-528(2004). Eichmueller S., Proc. Natl. Acad. Sci. U.S.A. 98:629-634(2001).