

## LMO3 Antibody [#C11382]

<b>Catalog Number:</b>	C11382
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q8TAP4
<b>Other Names:</b>	LIM domain only protein 3, Neuronal-specific transcription factor DAT1, Rhombotin-3, LMO3, RBTN3, RBTNL2, RHOM3
<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 LMO3.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	LMO3 antibody detects endogenous levels of total LMO3 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000   ELISA: 1:10000
<b>References:</b>	Ota T., Nat. Genet. 36:40-45(2004). The MGC Project Team, Genome Res. 14:2121-2127(2004). Foroni L., J. Mol. Biol. 226:747-761(1992).