

## TGF $\beta$ Receptor III Antibody [#B1128]

<b>Catalog Number:</b>	B1128
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q03167
<b>Other Names:</b>	TGBR3, TGF-beta receptor type III, TGFR-3, betaglycan, transforming growth factor beta receptor III
<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 TGF $\beta$ Receptor III.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	TGF $\beta$ Receptor III antibody detects endogenous levels of total SEK1/MKK4/JNKK1 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 ELISA: 1:5000
<b>References:</b>	SE Jacobsen, J. Immunol., Nov 1993; 151: 4534. T Parekh, J. Immunol., Mar 1994; 152: 2456. RW Pelton, Development, Oct 1990; 110: 609.