

## TNXB Antibody [#C19096]

<b>Catalog Number:</b>	C19096
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P22105
<b>Other Names:</b>	Tenascin-X; TN-X; Hexabrachion-like protein; TNXB; HXBL, TNX, TNXB1, TNXB2, XB
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), 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 TNXB.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	TNXB antibody detects endogenous levels of total TNXB protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 85%, Positives = 92%)
<b>Applications:</b>	IHC: 1:50~1:100 ELISA: 1:10000
<b>References:</b>	Xie T., Genome Res. 13:2621-2636(2003). Bristow J., J. Cell Biol. 122:265-278(1993). Tee M.K., Genomics 28:171-178(1995).