

## Trk A (Ab-496) Antibody [#B0589]

<b>Catalog Number:</b>	B0589
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P04629
<b>Other Names:</b>	EC 2.7.10.1, High affinity nerve growth factor receptor precursor, NTRK1, Slow nerve growth factor receptor, TRK, TRK1 transforming tyrosine kinase protein, TRKA, Trk-A, kinase TrkA, p140-TrkA
<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 non-phosphopeptide derived from human Trk A around the phosphorylation site of tyrosine 496 (P-Q-Y <sup>P</sup> -F-S).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Trk A (Ab-496) antibody detects endogenous levels of total Trk A protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	E Lavenius. Cell Growth Differ., Jun 1995; 6: 727. F Miralles. J. Endocrinol., Mar 1998; 156: 431 - 439. Tim R. Bilderback. J. Biol. Chem., Jan 1999; 274: 257. Darren Esposito. J. Biol. Chem., Aug 2001; 276: 32687 - 32695.