

Trk A Antibody [#B0034]

Catalog Number:	B0034
Concentration:	1mg/ml
Swiss-Prot No.:	P04629
Other Names:	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
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized peptide derived from Internal of human Trk A.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Trk A antibody detects endogenous levels of total Trk A protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:100~1:500 ELISA: 1:10000
References:	Martin-Zanca D., Mol. Cell. Biol. 9:24-33(1989). Shelton D.L., J. Neurosci. 15:477-491(1995). Indo Y., Jpn. J. Hum. Genet. 42:343-351(1997).