

Trk A Antibody [#B8070]

Catalog Number:	B8070
Concentration:	1mg/ml
Swiss-Prot No.:	P04629
Other Names:	EC 2.7.10.1; Neurotrophic tyrosine kinase receptor type 1; TRK1-transforming tyrosine kinase protein; p140-TrkA; Trk-A
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%); Mouse (Identities = 92%, Positives = 100%); Rat (Identities = 92%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:5000
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).