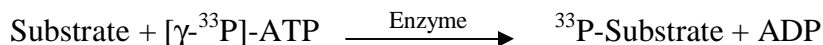


TrkB/NTRK2

(BDNF/NT-3 growth factors receptor [Precursor], Neurotrophic tyrosine kinase receptor type 2, TrkB tyrosine kinase, GP145-TrkB, NTRK2)

- **CAT#:** TrkB
- **Enzyme:** Human TrkB
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	31.5
PP2	173,000
PDGFR I-III	11,200
D-64406	2,500

- **Scientific Information:** Q16620 (<http://www.expasy.org/uniprot/Q16620>)

Reference Compound IC50 for TrkB

