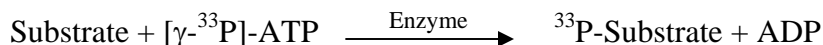


TRKB

(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), 0.2 mg/ml
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	<0.1
PP2	2,600
AG1478	15,000
PDGFR I-III	>20,000
D-64406	880

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

Reference compound IC50 for TRKB

