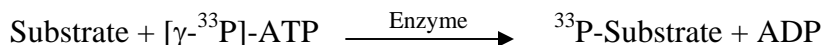


TRKC/NTRK3

(NT-3 growth factor receptor [Precursor], Neurotrophic tyrosine kinase receptor type 3, TrkC tyrosine kinase, GP145-TrkC, NTRK3)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	< 0.1
PP2	1,170
AG1478	1,700
PDGFR I-III	2,400
D-64406	129

- **Scientific Information:** Q16288 (<http://www.expasy.org/uniprot/Q16288>)

Reference compound IC50 for TRKC

