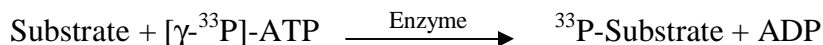


## TrkA/NTRK1

(High affinity nerve growth factor receptor [Precursor], Neurotrophic tyrosine kinase receptor type 1, TRK1 transforming tyrosine kinase protein, p140-TrkA, NTRK1)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.6
PP2	3,200
AG1478	6,200
PDGFR I-III	212
D-64406	328

- **Scientific Information:** P04629 (<http://www.expasy.org/uniprot/P04629>)

**Reference compound IC50 for TrkA**

