

TNK1

(Non-receptor tyrosine-protein kinase TNK1, CD38 negative kinase 1)

- **CAT#:** TNK1
- **Enzyme:** Human TNK1
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.4
PP2	20,000
AG1478	420
PDGFR I-III	8,400
D-64406	4,000

- **Scientific Information:** Q13470 (<http://www.expasy.org/uniprot/Q13470>)

Reference compound IC50 for TNK1

