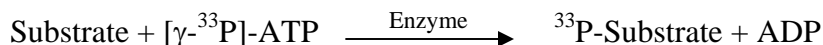


TIE2

(Angiopoietin-1 receptor [Precursor], Tyrosine-protein kinase receptor TIE-2, hTIE2, Tyrosine-protein kinase receptor TEK, p140 TEK, Tunica interna endothelial cell kinase)

- **CAT#:** TIE2
- **Enzyme:** Human TIE2
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	10
PP2	3,100
AG1478	5,300
D-64406	7,800

- **Scientific Information:** Q02763 (<http://www.expasy.org/uniprot/Q02763>)

Reference compound IC50 for TIE2

