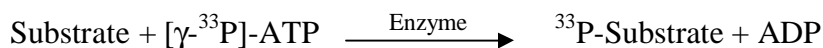


PYK2/FAK2/PTK2B

(Protein tyrosine kinase 2 beta, Focal adhesion kinase 2, FADK 2, Proline-rich tyrosine kinase 2, Cell adhesion kinase beta, CAK beta, Calcium-dependent tyrosine kinase, CADTK, Related adhesion focal tyrosine kinase, RAFTK, FAK2, PTK2B)

- **CAT#:** PYK2
- **Enzyme:** Human PYK2
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	2.7
PP2	10,700
AG1478	15,000
D-64406	29,000

- **Scientific Information:** Q14289 (<http://www.expasy.org/uniprot/Q14289>)

Reference compound IC50 for PYK2

