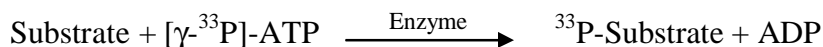


## FAK

(Focal adhesion kinase 1, FADK 1, pp125FAK, Protein-tyrosine kinase 2, PTK2, FAK1)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	33
PP2	895,000
AG1478	78,700
PDGFR I-III	45,000
D-64406	4,700

- **Scientific Information:** Q05397 (<http://www.expasy.org/uniprot/Q05397>)

**Reference Compound IC50 for FAK**

