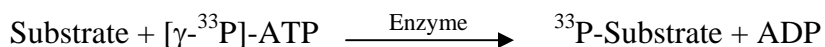


FER

(Proto-oncogene tyrosine-protein kinase FER, p94-FER, c-FER, Tyrosine kinase 3, TYK3)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.2

- **Scientific Information:** P16591 (<http://www.expasy.org/uniprot/P16591>)

Reference compound IC50 for FER

