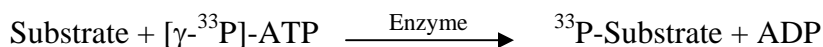


FYN

(Proto-oncogene tyrosine-protein kinase Fyn, p59-Fyn, Protooncogene Syn, SLK)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.9
PP2	20
AG1478	2,300
D-64406	500

- **Scientific Information:** P06241 (<http://www.expasy.org/uniprot/P06241>)

Reference compound IC50 for FYN

