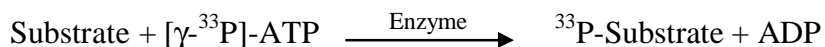


LYN

(Tyrosine-protein kinase Lyn)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.4
PP2	4.5
AG1478	62.8
PDGFR I-III	39,000
D-64406	470

- **Scientific Information:** P07948 (<http://www.expasy.org/uniprot/P07948>)

