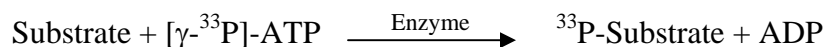


## LYN

(Tyrosine-protein kinase Lyn)

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

- **Reaction:**



- **Reference Compound IC50s:**

| Compounds     | IC50 (nM) |
|---------------|-----------|
| Staurosporine | 0.8       |
| PP2           | 46        |
| AG1478        | 260       |
| D-64406       | 630       |

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

### Reference compound IC50 for LYN

