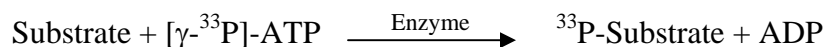


SYK

(Spleen tyrosine kinase, Tyrosine-protein kinase SYK, KSYK)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.3
AG1478	17,000
D-64406	10,000

- **Scientific Information:** P43405 (<http://www.expasy.org/uniprot/P43405>)

Reference compound IC50 for SYK

