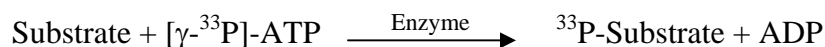


TYK1

(Leukocyte tyrosine kinase receptor [Precursor], Protein tyrosine kinase 1, LTK)

- **CAT#:** TYK1
- **Enzyme:** Human TYK1
- **Substrate:** Peptide substrate, [EAIYAAPFAKKK], 20 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	600
PP2	4,500
D-64406	18,100

- **Scientific Information:** P42680 (<http://www.expasy.org/uniprot/P42680>)

