

## TYK1

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

- **CAT#:** TYK1
- **Enzyme:** Human TYK1
- **Substrate:** Peptide substrate, [EAIYAAPFAKKK], 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	6.6
PP2	28,300
AG1478	1,100
D-64406	62,600

- **Scientific Information:** P29376 (<http://www.expasy.org/uniprot/P29376>)

Reference compound IC50 for TYK1/LTK

