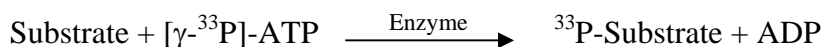


TYK2

(Non-receptor tyrosine-protein kinase TYK2)

- **CAT#:** TYK2
- **Enzyme:** Human TYK2
- **Substrate:** Peptide substrate, [KKSARGDYMTMQIG], 20 μ M
ATP 10 μ M
- **Reaction:**

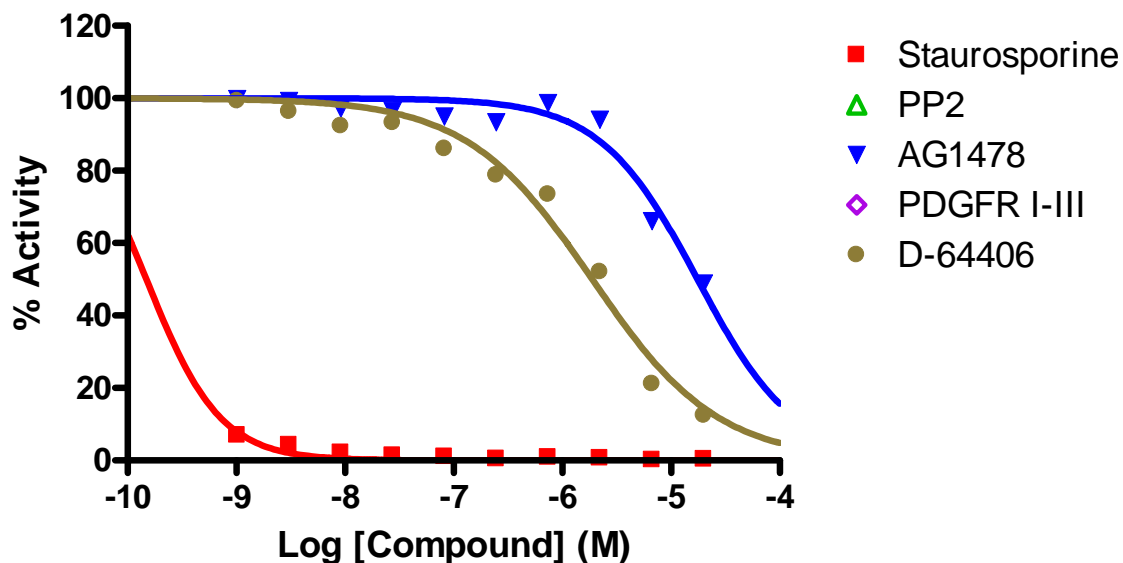


- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	< 0.1
AG1478	17,500
D-64406	1,900

- **Scientific Information:** P29597 (<http://www.expasy.org/uniprot/P29597>)

Reference compound IC50 for TYK2



Service performed
using kinases
under license from  **invitrogen**[™]
part of *life* technologies[™]