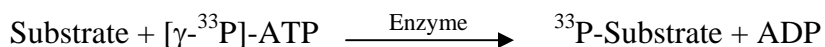


LCK

(Proto-oncogene tyrosine-protein kinase LCK, p56-LCK)

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

- **Reaction:**

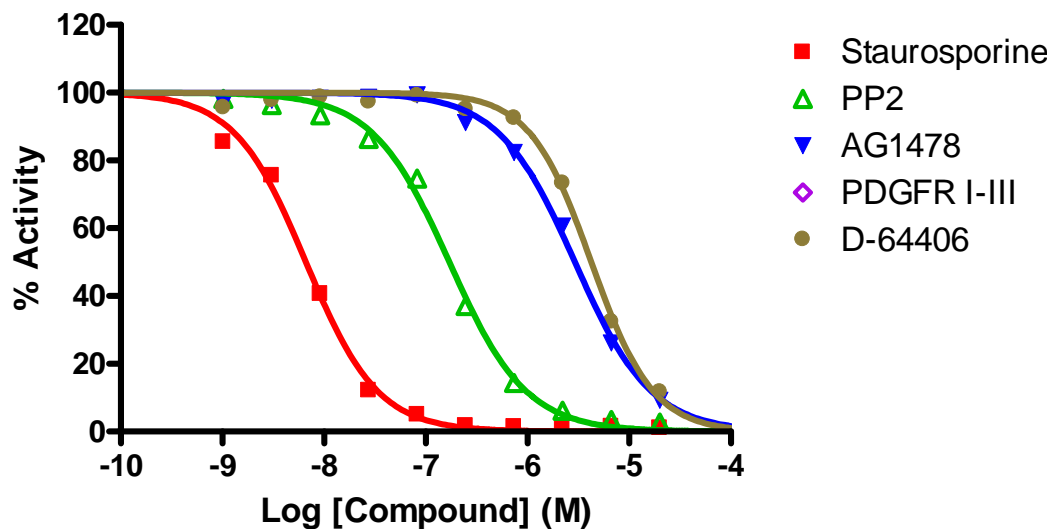


- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	6.6
PP2	170
AG1478	2,900
D-64406	4,300

- **Scientific Information:** P06239 (<http://www.expasy.org/uniprot/P06239>)

Reference compound IC50 for LCK



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