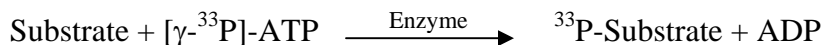


YES/YES1

(Proto-oncogene tyrosine-protein kinase Yes, p61-Yes, c-Yes)

- **CAT#:** YES/YES1
- **Enzyme:** Human YES/YES1
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml
ATP 10 μ M

- **Reaction:**

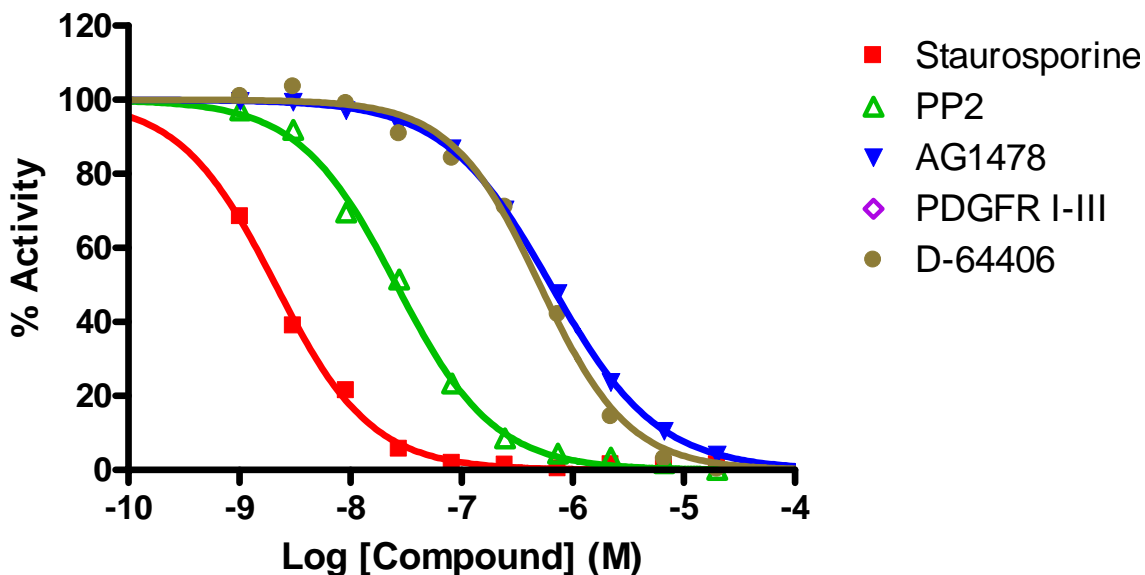


- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	2.1
PP2	26
AG1478	640
D-64406	510

- **Scientific Information:** P07947 (<http://www.expasy.org/uniprot/P07947>)

Reference compound IC50 for YES/YES1



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