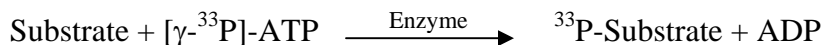


c-SRC

(Proto-oncogene tyrosine-protein kinase Src, pp60c-src)

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

- **Reaction:**

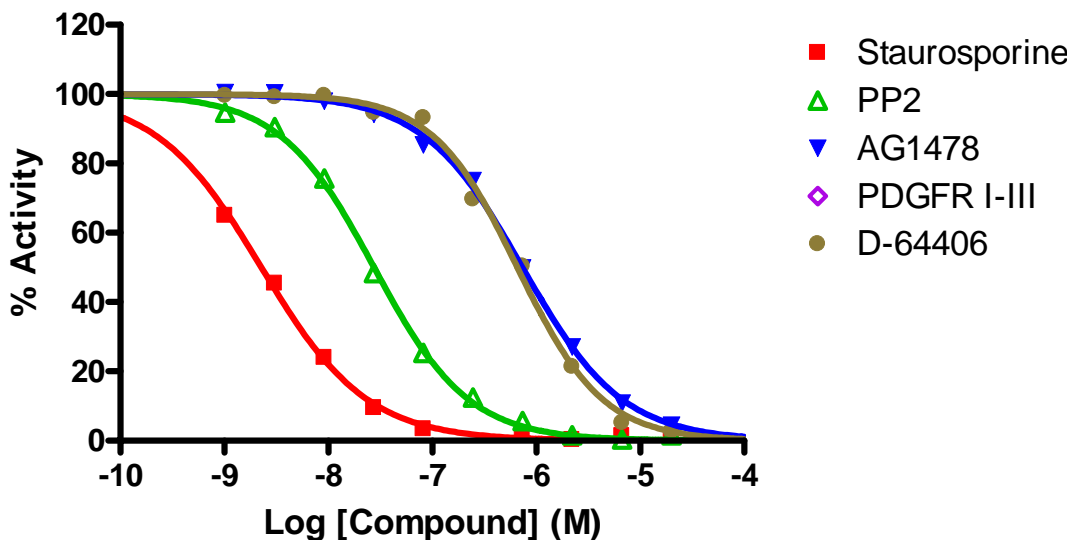


- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	2.3
PP2	28
AG1478	740
D-64406	660

- **Scientific Information:** P12931 (<http://www.expasy.org/uniprot/P12931>)

Reference compound IC50 for c-Src



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