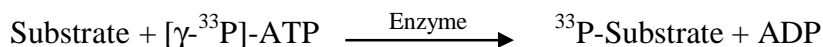


c-SRC

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

- **CAT#:** c-SRC
- **Enzyme:** Human c-SRC
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 1 μ M
- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	6.8
PP2	16.9
AG1478	590
D-64406	560

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

