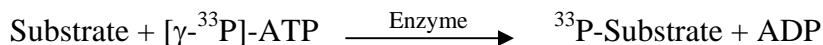


## c-SRC (T341M)

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

- **CAT#:** c-SRC (T341M)
- **Enzyme:** Human c-SRC (T341M)
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml  
ATP 1  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.24
D-64406	2,600

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

### Reference Compound IC50 for c-SRC (T341M)

