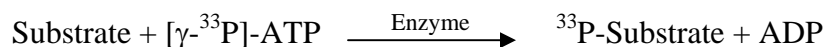


## CSK

(Tyrosine-protein kinase CSK, C-SRC kinase)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	14.6
PP2	142
AG1478	3,200
D-64406	9,500

- **Scientific Information:** P41240 (<http://www.expasy.org/uniprot/P41240>)

Reference compound IC50 for CSK

