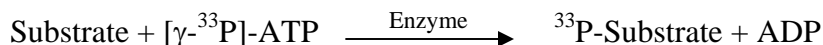


CTK

(Megakaryocyte-associated tyrosine-protein kinase, Tyrosine-protein kinase CTK, Protein kinase HYL, Hematopoietic consensus tyrosine-lacking kinase, MATK, HYL)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	770

- **Scientific Information:** P42679 (<http://www.expasy.org/uniprot/P42679>)

Reference compound IC50 for CTK/MATK/HYL

