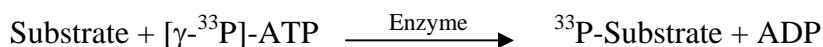


HCK

(Tyrosine-protein kinase HCK, p59-HCK/p60-HCK, Hemopoietic cell kinase)

- **CAT#:** HCK
- **Enzyme:** Human HCK
- **Substrate:** Peptide substrate, [KVEKIGEGTYGVVYK], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.7
PP2	46
AG1478	540
D-64406	1,400

- **Scientific Information:** P08631 (<http://www.expasy.org/uniprot/P08631>)

Reference compound IC50 for HCK

