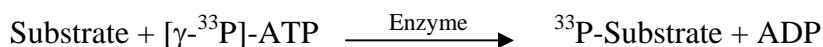


CHK2

(Serine/threonine-protein kinase Chk2, Cds1, CHEK2)

- **CAT#:** CHK2
- **Enzyme:** Human CHK2
- **Substrate:** Peptide substrate, [KKKVSRSGLYRSPSPENLNRPR], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	5.2
Ro 31-8220	690
GW 5074	1,200
H-89	2,500
Rottlerin	3,300

- **Scientific Information:** O96017 (<http://www.expasy.org/uniprot/O96017>)

Reference compound IC50 for CHK2

