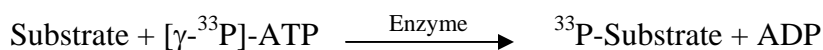


CDK9/Cyclin K

(Cell division protein kinase 9, Cyclin-dependent kinase 9, Serine/threonine-protein kinase PITALRE, C-2K, Cell division cycle 2-like protein kinase 4, CDC2L4
Cyclin K: Cyclin-K, CCNK, CPR4)

- **CAT#:** CDK9/Cyclin K
- **Enzyme:** Human CDK9/Cyclin K
- **Substrate:** Peptide substrate, [KTFCGTPEYLAPEVRREPRILSEEEQEMFRDFDYIADWC], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	11
Ro 31-8220	8.8
GW 5074	5,100
H-89	10,800
Rottlerin	17,700

- **Scientific Information:** CDK9: P50750 (<http://www.expasy.org/uniprot/P50750>)
Cyclin K: O75909 (<http://www.expasy.org/uniprot/O75909>)

Reference Compound IC50 for CDK 9/cyclin K

