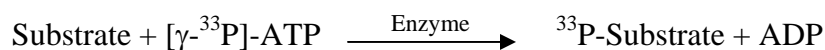


## CDK1/Cyclin E

(CDK1: Cyclin-dependent kinase 1, p34 protein kinase, CDC2,  
Cell division control protein 2 homolog,  
Cyclin E: G1/S-specific cyclin-E1, CCNE1)

- **CAT#:** CDK1/Cyclin E
- **Enzyme:** Human CDK1/Cyclin E
- **Substrate:** RB protein, 1  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	3.8
Ro 31-8220	13.2
GW 5074	10,000
H-89	20,000
Rottlerin	>50,000

- **Scientific Information:** CDK1: P06493 (<http://www.expasy.org/uniprot/P06493>)  
Cyclin E: P24864 (<http://www.expasy.org/uniprot/P24864>)

### Reference compound IC50 for CDK1/cyclin E

