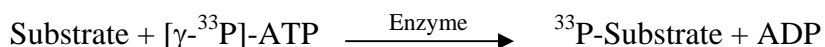


## CDK2/Cyclin A1

(CDK2: Cyclin-dependent kinase 2, p33 protein kinase, Cell division protein kinase 2,  
Cyclin A1: CCNA1)

- **CAT#:** CDK2/Cyclin A1
- **Enzyme:** Human CDK2/Cyclin A1
- **Substrate:** Histone H1, 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

| Compounds     | IC50 (nM) |
|---------------|-----------|
| Staurosporine | 1.2       |
| Ro 31-8220    | 4.6       |
| GW 5074       | >50,000   |
| H-89          | 11,000    |
| Rottlerin     | >50,000   |

- **Scientific Information:** CDK2: P24941 (<http://www.expasy.org/uniprot/P24941>)  
Cyclin A1: P78396 (<http://www.expasy.org/uniprot/P78396>)

### Reference compound IC50 for CDK2/cyclin A1

