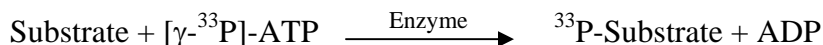


CDK2/Cyclin A2

(CDK2: Cyclin-dependent kinase 2, Cell division protein kinase 2, p33 protein kinase,
Cyclin A2: CCNA2)

- **CAT#:** CDK2/Cyclin A2
- **Enzyme:** Human CDK2/Cyclin A2
- **Substrate:** Histone H1, 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.3
Ro 31-8220	3.7
GW 5074	5,700
H-89	12,000

- **Scientific Information:** CDK2: P24941 (<http://www.expasy.org/uniprot/P24941>)
Cyclin A2: P20248 (<http://www.expasy.org/uniprot/P20248>)

Reference compound IC50 for CDK2/cyclin A

