

CDK1/Cyclin A

(CDK1: Cyclin-dependent kinase 1, p34 protein kinase, CDC2,
Cell division control protein 2 homolog
Cyclin A: Cyclin-A2, Cyclin-A, CCNA2, CCN1, CCNA)

- **CAT#:** CDK1/cyclin A
- **Enzyme:** Human CDK1/cyclin A
- **Substrate:** Histone H1, 20 μ M
ATP 1 μ M
- **Reaction:**

$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} \text{}^{33}\text{P-Substrate} + \text{ADP}$$
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	6.6
Ro 31-8220	58.4
H-89	42,000

- **Scientific Information:** CDK1: P06493 (<http://www.expasy.org/uniprot/P06493>)
Cyclin A: P20248 (<http://www.expasy.org/uniprot/P20248>)

Reference Compound IC50 for CDK1/Cyclin A

