

CDK3/Cyclin E

(CDK3: Cell division protein kinase 3, Cyclin-dependent kinase 3,
Cyclin E: G1/S-specific cyclin-E1, CCNE1)

- **CAT#:** CDK3/Cyclin E
- **Enzyme:** Human CDK3/Cyclin E
- **Substrate:** Histone H1, 5 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	1.2
Ro 31-8220	3.3
GW5074	22,000
H-89	11,000

- **Scientific Information:** CDK3: Q00526 (<http://www.expasy.org/uniprot/Q00526>)
Cyclin E: P24864 (<http://www.expasy.org/uniprot/P24864>)

Reference compound IC50 for CDK3/Cyclin E

