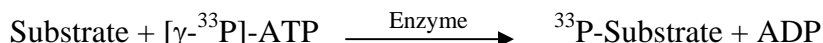


CDK4/Cyclin D3

(CDK4: Cell division protein kinase 4, Cyclin-dependent kinase 4, PSK-J3
Cyclin D3: G1/S-specific cyclin-D3, CCND3)

- **CAT#:** CDK4/Cyclin D3
- **Enzyme:** Human CDK4/Cyclin D3
- **Substrate:** RB-CTF, 1 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	83
Ro 31-8220	25.8
GW 5074	21,500
H-89	213,000
Rottlerin	227,000

- **Scientific Information:** CDK4: P11802 (<http://www.expasy.org/uniprot/P11802>)
Cyclin D3: P30281 (<http://www.expasy.org/uniprot/P30281>)

Reference Compound IC50 for CDK4/Cylin D3

