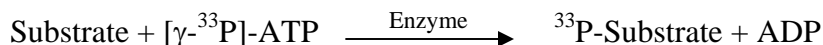


## PLK2

(Serine/threonine-protein kinase PLK2, Polo-like kinase 1, PLK-2,  
Serine/threonine-protein kinase SNK, Serum-inducible kinase, SNK)

- **CAT#:** PLK2
- **Enzyme:** Human PLK2
- **Substrate:** Casein, 1 mg/ml  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	740
Ro 31-8220	49,000
GW 5074	19,000

- **Scientific Information:** Q9NYY3 (<http://www.expasy.org/uniprot/Q9NYY3>)

### Reference compound IC50 for PLK2

