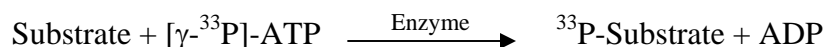


## PLK4/SAK

(Serine/threonine-protein kinase PLK4, Polo-like kinase 4,  
Serine/threonine-protein kinase 18, Serine/threonine-protein kinase Sak, STK18)

- **CAT#:** PLK4/SAK
- **Enzyme:** Human PLK4/SAK
- **Substrate:** Casein, 2 mg/ml  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	4.2
Ro 31-8220	15,000
GW5074	4,700
H-89	>50,000
Rottlerin	>50,000

- **Scientific Information:** O00444 (<http://www.expasy.org/uniprot/O00444>)

### Reference compound IC50 for PLK4/SAK

