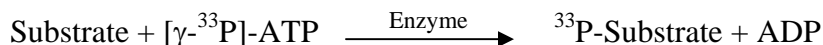


## SLK/STK2

(STE20-like serine/threonine-protein kinase, STE20-related serine/threonine-protein kinase, STE20-related kinase, Serine/threonine-protein kinase 2, CTCL tumor antigen se20-9, KIAA0204)

- **CAT#:** SLK
- **Enzyme:** Human SLK
- **Substrate:** Histone H3, 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	3.9
Ro 31-8220	38.9
GW5074	18,000
H-89	42,000

- **Scientific Information:** Q9H2G2 (<http://www.expasy.org/uniprot/Q9H2G2>)

### Reference compound IC50 for SLK/STK2

