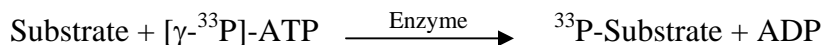


CLK4

(Dual specificity protein kinase CLK4, CDC-like kinase 4)

- **CAT#:** CLK4
- **Enzyme:** Human CLK4
- **Substrate:** MBP (Myelin Basic Protein), 10 μ M
ATP 10 μ M

- **Reaction:**

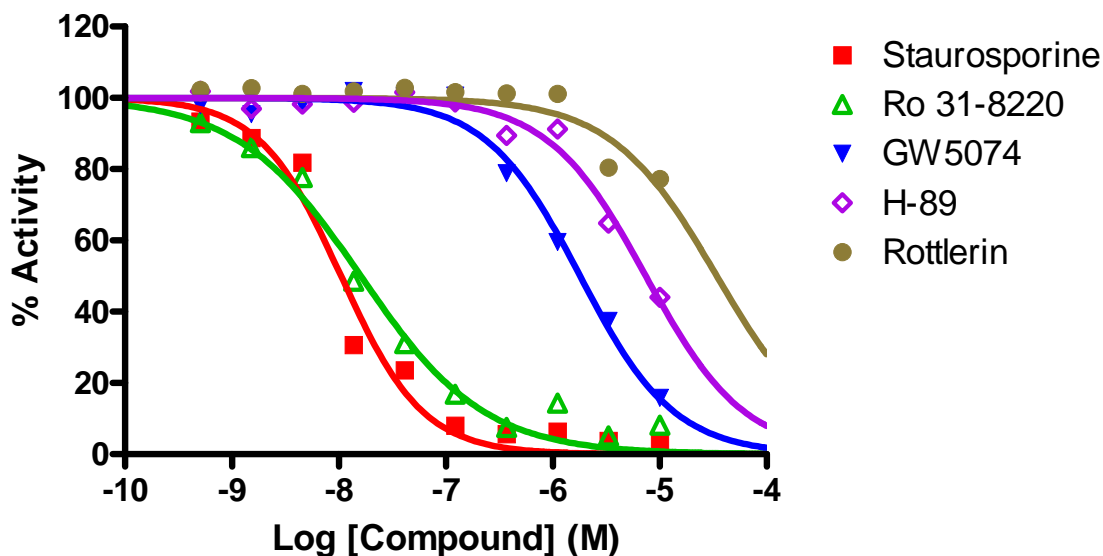


- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	10.5
Ro 31-8220	16
GW 5074	1,800
H-89	7,400
Rottlerin	34,000

- **Scientific Information:** Q9HAZ1 (<http://www.expasy.org/uniprot/Q9HAZ1>)

Reference compound IC50 for CLK4



Service performed
using kinases
under license from  **invitrogen**[™]
part of *life* technologies[™]