

SIK2

(Serine/threonine-protein kinase SNF1-like kinase 2, Qin-induced kinase
Salt-inducible protein kinase 2, SNF1LK2, KIAA0781, QIK, SN1L2)

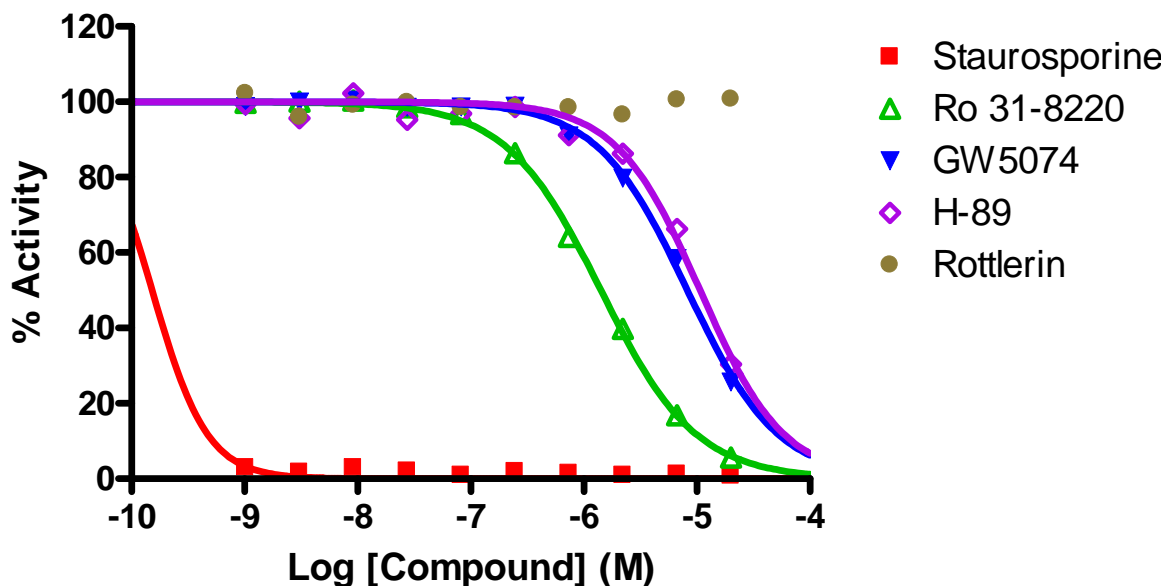
- **CAT#:** SIK2
- **Enzyme:** Human SIK2
- **Substrate:** Peptide substrate, [AMARAASAAALARRR], 20 μ M
ATP 10 μ M
- **Reaction:**

$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} \text{}^{33}\text{P-Substrate} + \text{ADP}$$
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	<0.1
Ro 31-8220	1,400
GW5074	8,300
H-89	11,000
Rottlerin	>20,000

- **Scientific Information:** Q9H0K1 (<http://www.expasy.org/uniprot/Q9H0K1>)

Reference compound IC50 for SIK2



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