

SIK1

(Serine/threonine-protein kinase SIK1, Salt-inducible protein kinase 1,
Serine/threonine-protein kinase SNF1-like kinase 1, SNF1LK, SIK)

- **CAT#:** SIK1
- **Enzyme:** Human SIK1
- **Substrate:** Peptide substrate, [AMARAASAAALARRR], 20 μ M
ATP 10 μ M
- **Reaction:**

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

Compounds	IC50 (nM)
Staurosporine	1.2
Ro 31-8220	3,000
GW5074	3,600
H-89	29,000
Rottlerin	>30,000

- **Scientific Information:** P57059 (<http://www.expasy.org/uniprot/P57059>)

Reference compound IC50 for SIK1

