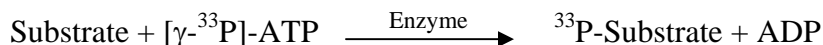


## SGK3

(Serine/threonine-protein kinase Sgk3, Serum/glucocorticoid-regulated kinase 3  
Serum/glucocorticoid-regulated kinase-like, SGK2, SGKL)

- **CAT#:** SGK3
- **Enzyme:** Human SGK3
- **Substrate:** Peptide substrate, [GRPRTSSFAEG], 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	236
Ro 31-8220	745
Rottlerin	2,300

- **Scientific Information:** Q96BR1 (<http://www.expasy.org/uniprot/Q96BR1>)

### Reference compound IC50 for SGK3/SGKL

