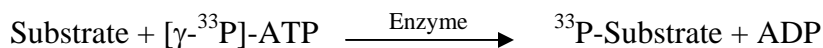


MSK2

(Ribosomal protein S6 kinase alpha-4, Nuclear mitogen- and stress-activated protein kinase 2, 90 kDa ribosomal protein S6 kinase 4, Ribosomal protein kinase B, RSKB, RPS6KA4, KS6A4)

- **CAT#:** MSK2
- **Enzyme:** Human MSK2
- **Substrate:** Peptide substrate, [GRPRTSSFAEG], 20 μ M
ATP 10 μ M

- **Reaction:**

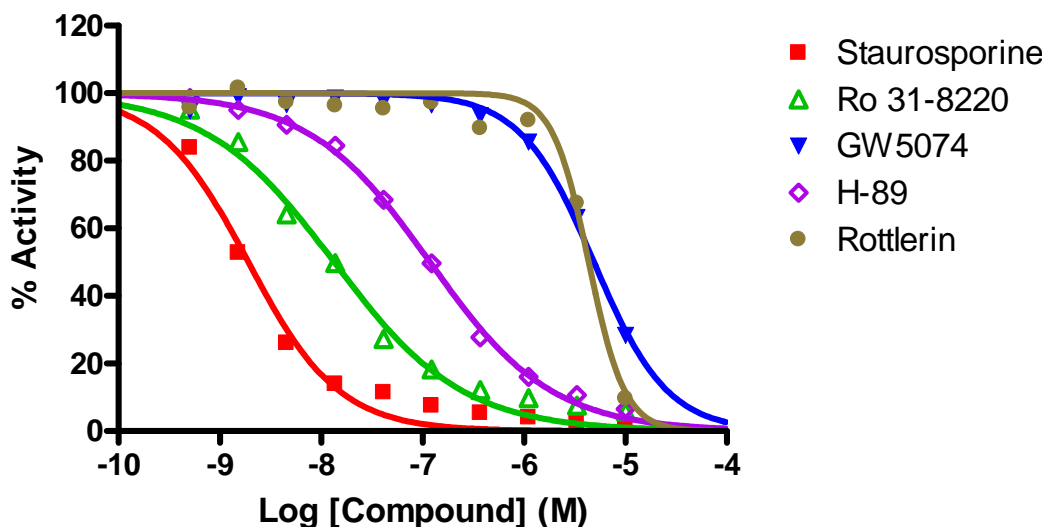


- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.9
Ro 31-8220	14
GW 5074	4,900
H-89	120
Rottlerin	4,400

- **Scientific Information:** O75676 (<http://www.expasy.org/uniprot/O75676>)

Reference Compound IC50 for MSK2



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