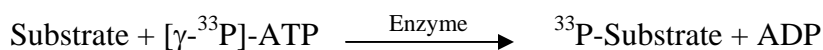


## p70S6Kb/RPS6KB2

(Ribosomal protein S6 kinase beta-2, S6K-beta 2, S6K2, 70 kDa ribosomal protein S6 kinase 2, p70 ribosomal S6 kinase beta, p70 S6 kinase beta, p70 S6Kbeta, p70-S6KB, S6K-beta, p70-beta, S6 kinase-related kinase, SRK, Serine/threonine-protein kinase 14 beta, RPS6KB2, STK14B, KS6B2)

- **CAT#:** p70S6Kb/RPS6KB2
- **Enzyme:** Human p70S6Kb/RPS6KB2
- **Substrate:** Peptide substrate, [KKRNRTLTK], 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	4.9
Ro 31-8220	38
GW 5074	14,000
H-89	1,400
Rottlerin	12,000

- **Scientific Information:** Q9UBS0 (<http://www.expasy.org/uniprot/Q9UBS0>)

### Reference Compound IC50 for p70S6Kb/RPS6KB2

