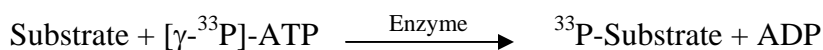


p70S6K/RPS6KB1

(Ribosomal protein S6 kinase beta-1, Ribosomal protein S6 kinase I, S6K, S6K1, 70 kDa ribosomal protein S6 kinase 1, p70 S6 kinase alpha, p70(S6K)-alpha, p70-S6K, P70S6K, p70-alpha, KS6B1, RPS6KB1)

- **CAT#:** p70S6K/RPS6KB1
- **Enzyme:** Human p70S6K/RPS6KB1
- **Substrate:** Peptide substrate, [KKRNRTLTK], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.9
Ro 31-8220	2.8
GW 5074	5,100
H-89	140
Rottlerin	37,000

- **Scientific Information:** P23443 (<http://www.expasy.org/uniprot/P23443>)

Reference compound IC50 for p70S6K/RPS6KB1

