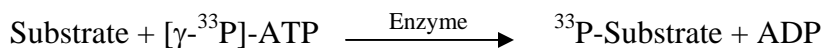


mTOR

(FKBP12-rapamycin complex-associated protein, Rapamycin target protein, RAPT1, FK506-binding protein 12-rapamycin complex-associated protein 1, Mammalian target of rapamycin, FRAP, FRAP1, FRAP2)

- **CAT#:** mTOR
- **Enzyme:** Human mTOR
- **Substrate:** p70S6K, 0.5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	91
Ro 31-8220	390
GW 5074	1,800
H-89	131,000

- **Scientific Information:** P42345 (<http://www.expasy.org/uniprot/P42345>)

Reference Compound IC50 for mTOR

