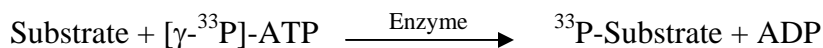


## mTOR

(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:** 4EBP1 1  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**

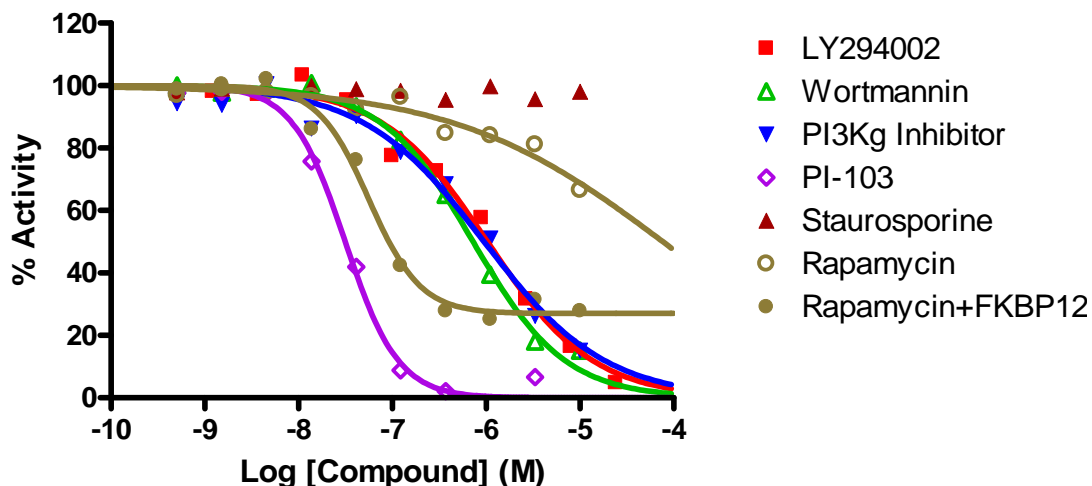


- **Reference Compound IC50s:**

Compounds	IC50 (nM)
LY294002	1,030
Wortmannin	727
PI3K $\gamma$ Inhibitor	952
PI-103	31.5
Rapamycin	77,000
Rapamycin + FKBP12	56

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

### Reference Compound IC50 for mTOR



Service performed using kinases under license from  **invitrogen**<sup>™</sup>  
part of *life* technologies<sup>™</sup>