

TTBK2

(Tau-tubulin kinase 2, KIAA0847)

- **CAT#:** TTBK2
- **Enzyme:** Human TTBK2
- **Substrate:** MBP (Myelin Basic Protein), 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	77,000
Ro 31-8220	>100,000
GW 5074	>100,000
H-89	>100,000
Rottlerin	>100,000

- **Scientific Information:** Q6IQ55 (<http://www.expasy.org/uniprot/Q6IQ55>)

Reference compound IC50 for TTBK2

