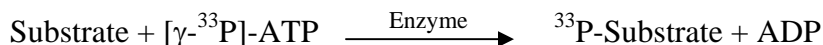


TAOK1

(Serine/threonine-protein kinase TAO1, Thousand and one amino acid protein 1, STE20-like kinase PSK2, Kinase from chicken homolog B, hKFC-B, KIAA1361, MAP3K16)

- **CAT#:** TAOK1
- **Enzyme:** Human TAOK1
- **Substrate:** MBP (Myelin Basic Protein), 10 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.4
Ro 31-8220	186
GW 5074	2,100
H-89	13,500
Rottlerin	28,500

- **Scientific Information:** Q7L7X3 (<http://www.expasy.org/uniprot/Q7L7X3>)

Reference Compound IC50 for TAOK1

