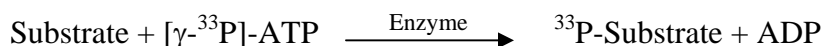


CAMKK2

(Calcium/calmodulin-dependent protein kinase kinase 2, Calcium/calmodulin-dependent protein kinase, kinase beta, CaM-kinase kinase beta, CaM-KK beta, CaMKK beta, KKCC2, CAMKKB, KIAA0787)

- **CAT#:** CAMKK2
- **Enzyme:** Human CAMKK2
- **Substrate:** MBP (Myelin Basic Protein), 10 μ M
Ca²⁺-Calmodulin, 1 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	17.4
Ro 31-8220	147
GW 5074	1,000

- **Scientific Information:** Q96RR4 (<http://www.expasy.org/uniprot/Q96RR4>)

Reference Compound IC50 for CAMKK2

