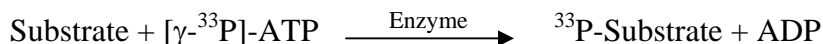


CAMK2a

(Calcium/calmodulin-dependent protein kinase type II alpha chain, CaM-kinase II appha chain, CaM kinase II alpha subunit, CaMK-II subunit alpha, CAMK2A, KCC2A)

- **CAT#:** CAMK2a
- **Enzyme:** Human CAMK2a
- **Substrate:** Peptide substrate, [KKLNRTLSEFAEPG], 20 μ M
Ca²⁺-Calmodulin, 1 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	< 0.1
Ro 31-8220	6,800
GW 5074	21,600

- **Scientific Information:** Q9UQM7 (<http://www.expasy.org/uniprot/Q9UQM7>)

Reference Compound IC50 for CAMK2 α

