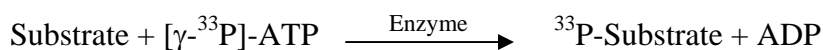


CAMKII γ

(Calcium/calmodulin-dependent protein kinase type II gamma chain, CaM-kinase II gamma chain
CaM kinase II gamma subunit, CaMK-II subunit gamma, CAMK2G, KCC2G)

- **CAT#:** CAMKII γ
- **Enzyme:** Human CAMKII γ
- **Substrate:** Peptide substrate, [KKLNRTLSFAEPG], 5 μ M
Ca²⁺-Calmodulin, 1 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	< 0.5
Ro 31-8220	544
GW 5074	3,600

- **Scientific Information:** Q13555 (<http://www.expasy.org/uniprot/Q13555>)

Reference Compound IC50 for CAMKII γ

