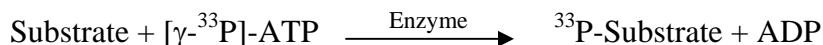


PHK γ 1

(Phosphorylase b kinase gamma catalytic chain, skeletal muscle isoform,
Phosphorylase kinase subunit gamma 1, PHKG1, PHKG)

- **CAT#:** PHK γ 1
- **Enzyme:** Human PHK γ 1
- **Substrate:** Peptide substrate, [KKLNRTLSFAEPG], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	3.3
Ro 31-8220	240
GW5074	11,000
H-89	12,400

- **Scientific Information:** Q16816 (<http://www.expasy.org/uniprot/Q16816>)

Reference compound IC50 for PHK γ 1

