

PHKg2

(Phosphorylase b kinase gamma catalytic chain, testis/liver isoform,
Phosphorylase kinase subunit gamma-2, PSK-C3, PHKG2)

- **CAT#:** PHKg2
- **Enzyme:** Human PHKg2
- **Substrate:** Peptide substrate, [KKLNRTLSEFAEPG], 20 μM
ATP 10 μM
- **Reaction:**

$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} \text{}^{33}\text{P}\text{-Substrate} + \text{ADP}$$
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.4
Ro 31-8220	22,000
GW 5074	9,400
H-89	8,900

- **Scientific Information:** P15735 (<http://www.expasy.org/uniprot/P15735>)

Reference compound IC50 for PHKg2

