

PDK1/PDPK1

(3-phosphoinositide-dependent protein kinase 1, hPDK1, PDK1)

- **CAT#:** PDK1/PDPK1
- **Enzyme:** Human PDK1/PDPK1
- **Substrate:** Peptide substrate, [KTFCGTPEYLAPEVRREPRILSEEEQEMFRDFDYIADWC], 20 μ M
ATP 10 μ M
- **Reaction:**

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

Compounds	IC50 (nM)
Staurosporine	13
Ro 31-8220	430
GW 5074	>20,000

- **Scientific Information:** O15530 (<http://www.expasy.org/uniprot/O15530>)

Reference compound IC50 for PDK1/PDPK1

