

## PDK1/PDHK1

(Pyruvate dehydrogenase [lipoamide]] kinase isozyme 1 mitochondrial,  
Pyruvate dehydrogenase kinase isoform 1)

- **CAT#:** PDK1/PDHK1
- **Enzyme:** Human PDK1/PDHK1
- **Substrate:** MBP (Myelin Basic Protein), 20  $\mu$ M  
ATP 10  $\mu$ 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	>100,000
Ro 31-8220	>100,000
GW 5074	9,000
H-89	>100,000
Rottlerin	>100,000

- **Scientific Information:** Q15118 (<http://www.expasy.org/uniprot/Q15118>)

### Reference compound IC50 for PDK1/PDHK1

