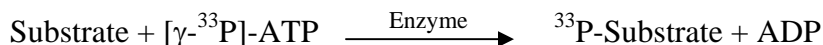


PDK3/PDHK3

(Pyruvate dehydrogenase [lipoamide]] kinase isozyme 3 mitochondrial,
Pyruvate dehydrogenase kinase isoform 3)

- **CAT#:** PDK3/PDHK3
- **Enzyme:** Human PDK3/PDHK3
- **Substrate:** MBP (Myelin Basic Protein), 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	>100,000
Ro 31-8220	>100,000
GW 5074	4,500
H-89	>100,000
Rottlerin	>100,000

- **Scientific Information:** Q15120 (<http://www.expasy.org/uniprot/Q15120>)

Reference compound IC50 for PDK3/PDHK3

