

## PDK4/PDKH4

([Pyruvate dehydrogenase [lipoamide]] kinase isozyme 4, mitochondrial,  
Pyruvate dehydrogenase kinase isoform 4)

- **CAT#:** PDK4/PDKH4
- **Enzyme:** Human PDK4/PDKH4
- **Substrate:** MBP (Myelin Basic Protein), 10  $\mu$ M  
ATP 10  $\mu$ M
- **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) |
|------------|-----------|
| Ro 31-8220 | >100,000  |
| GW 5074    | 5,000     |
| Rottlerin  | 47,600    |

- **Scientific Information:** Q16654 (<http://www.expasy.org/uniprot/Q16654>)

### Reference compound IC50 for PDK4/PDKH4

