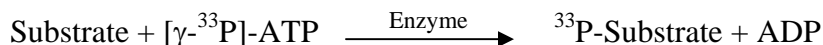


## EIF2AK3/PERK

(Eukaryotic translation initiation factor 2-alpha kinase 3, PRKR-like endoplasmic reticulum kinase, Pancreatic eIF2-alpha kinase, HsPEK, PEK, E2AK3)

- **CAT#:** EIF2AK3
- **Enzyme:** Human EIF2AK3
- **Substrate:** MBP (Myelin Basic Protein), 10  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	12,000
Ro 31-8220	27,300
GW 5074	17,700
Rottlerin	7,500

- **Scientific Information:** Q9NZJ5 (<http://www.expasy.org/uniprot/Q9NZJ5>)

### Reference compound IC50 for EIF2AK3/PERK

