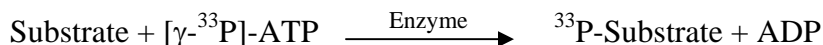


EIF2AK4/GCN2

(Eukaryotic translation initiation factor 2-alpha kinase 4, GCN2-like protein, KIAA1338, E2AK4)

- **CAT#:** EIF2AK4/GCN2
- **Enzyme:** Human EIF2AK4/GCN2
- **Substrate:** Peptide substrate, [GRSRSRSRSR], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	700
Ro 31-8220	>50,000
GW 5074	>50,000
H-89	>50,000
Rottlerin	>50,000

- **Scientific Information:** Q9P2K8 (<http://www.expasy.org/uniprot/Q9P2K8>)

Reference compound IC50 for EIF2AK4/GCN2

