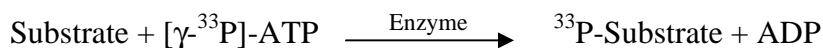


## AMPK (A2/B2/G2)

(Subunit A2: 5'-AMP-activated protein kinase catalytic subunit alpha-2, PRKAA2, AMPK2,  
Subunit B2: PRKAB2, AMPKb2, Subunit G2: PRKAG2, AMPKg2, H91620p)

- **CAT#:** AMPK (A2/B2/G2)
- **Enzyme:** Human AMPK (combination of A2/B2/G2 subunits)
- **Substrate:** Peptide substrate, [HMRSAMSGHLVKRR], 20  $\mu$ M  
AMP, 125  $\mu$ M  
ATP 1  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	<0.1
Ro 31-8220	100
GW 5074	2,800
H-89	990

- **Scientific Information:** Subunit A2: P54646 (<http://www.expasy.org/uniprot/P54646>)  
Subunit B2: O43741 (<http://www.expasy.org/uniprot/O43741>)  
Subunit G2: Q9UGJ0 (<http://www.expasy.org/uniprot/Q9UGJ0>)

### Reference compound IC50 for AMPK(A2/B2/G2)

