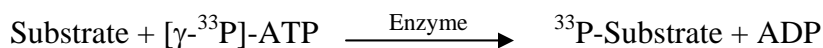


## AMPK (A2/B1/G1)

(Subunit A2: 5'-AMP-activated protein kinase catalytic subunit alpha-2, PRKAA2, AMPK2,  
Subunit B1: PRKAB1, AMPK, AMPKb, Subunit G1: PRKAG1, AMPKg)

- **CAT#:** AMPK (A2/B1/G1)
- **Enzyme:** Human AMPK (combination of A2/B1/G1 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	170
GW 5074	2,000
H-89	1,300
Rottlerin	50,300

- **Scientific Information:** Subunit A2: P54646 (<http://www.expasy.org/uniprot/P54646>)  
Subunit B1: Q9Y478 (<http://www.expasy.org/uniprot/Q9Y478>)  
Subunit G1: P54619 (<http://www.expasy.org/uniprot/P54619>)

### Reference Compound IC50 for AMPK (A2/B1/G1)

