

---

## AMPK $\alpha$ 2/ $\beta$ 1/ $\gamma$ 1(PRKAA2/B1/G1)

Product code 02-114

Full-length human AMPK $\alpha$ 2 [1-552(end) amino acids of accession number NP\_006243.2] was co-expressed as N-terminal GST-fusion protein (89 kDa) with GST-PRKAB1 [1-270(end) amino acids of accession number NP\_006244.2] and PRKAG1 [1-331(end) amino acids of accession number NP\_002724.1] using baculovirus expression system. GST-AMPK $\alpha$ 2/ $\beta$ 1/ $\gamma$ 1 was purified by using glutathione sepharose chromatography and activated with His-tagged CaMKK1. Activated GST-AMPK $\alpha$ 2/ $\beta$ 1/ $\gamma$ 1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay  
Substrate : SAMS peptide

ATP ( $\mu$ M) Km app / Bin : 100 / 100  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 0.79  
IC50 at 1 mM ATP (nM) : n.a.

---