

# AMPK $\alpha$ 1/ $\beta$ 1/ $\gamma$ 1(PRKAA1/B1/G1)

Product code: 02-113

Full-length human AMPK $\alpha$ 1 [1-550(end) amino acids of accession number NP\_006242.4] was co-expressed as N-terminal GST-fusion protein (90 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 $\alpha$ 1/ $\beta$ 1/ $\gamma$ 1 was purified by using glutathione sepharose chromatography and activated with His-tagged CaMKK1. Activated GST-AMPK $\alpha$ 1/ $\beta$ 1/ $\gamma$ 1 was purified by using glutathione sepharose chromatography.

Assay platform	: MSA
Substrate	: SAMS peptide
ATP ( $\mu$ M) Km app / Bin	: 130 / 150
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.41
IC50 at 1 mM ATP (nM)	: 0.87