

AMPK α 1/ β 1/ γ 1(PRKAA1/B1/G1)

Product code: 02-113

Truncated human AMPK α 1 [10-559(end) amino acids of accession number NP_006242.5] 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 α 1/ β 1/ γ 1 was purified by using glutathione sepharose chromatography and activated with His-tagged CaMKK1. Activated GST-AMPK α 1/ β 1/ γ 1 was purified by using glutathione sepharose chromatography.

Assay platform	: MSA
Substrate	: SAMS peptide
ATP (μ 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