
AMPK α 1/ β 1/ γ 1(PRKAA1/B1/G1)

Product code 02-113

Full-length human AMPK α 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 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 : Mobility Shift Assay
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
