

Product Information

BTN-AMPK α 1/ β 2/ γ 1(PRKAA1/B2/G1)

Product Number : **02-447-20N**

Product description

Full-length human AMPK α 1 [1-550(end) amino acids of accession number NP_006242.4] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (67 kDa) with GST-PRKAB2 [1-272(end) amino acids of accession number NP_005390.1] and PRKAG1 [1-331(end) amino acids of accession number NP_002724.1] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose and activated with His-tagged CaMKK1. Activated protein was purified by using DYKDDDDK tag antibody agarose.

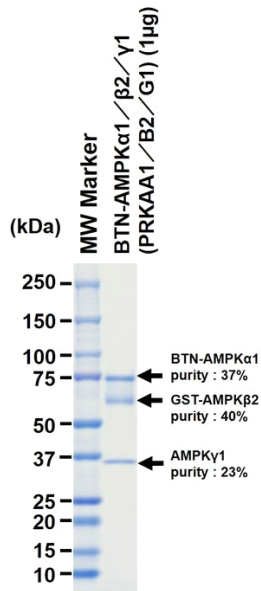
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
 1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

Store at -80C.
 Avoid repeating freeze-thaws.

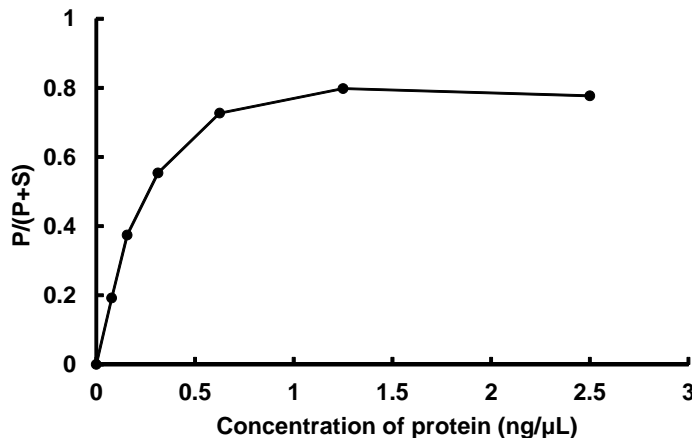
SDS-PAGE



Purity: 100 %

The purity was assessed by SDS-PAGE/CBB staining.

Activity data



The activity was measured by off-chip mobility shift assay(MSA). The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : SAMS peptide

ATP : 100 μ M