

Product Information

BTN-AMPK α 1/ β 1/ γ 1(PRKAA1/B1/G1)

Product Number : **02-413-20N**

Product description

Truncated human AMPK α 1 [10-559(end) amino acids of accession number NP_006242.5] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (67 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. 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.

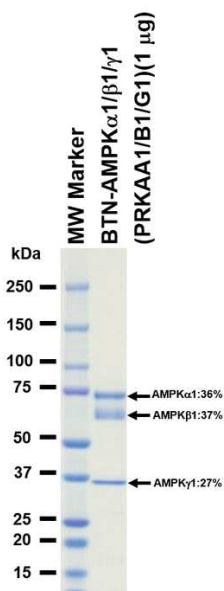
Storage buffer:

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.

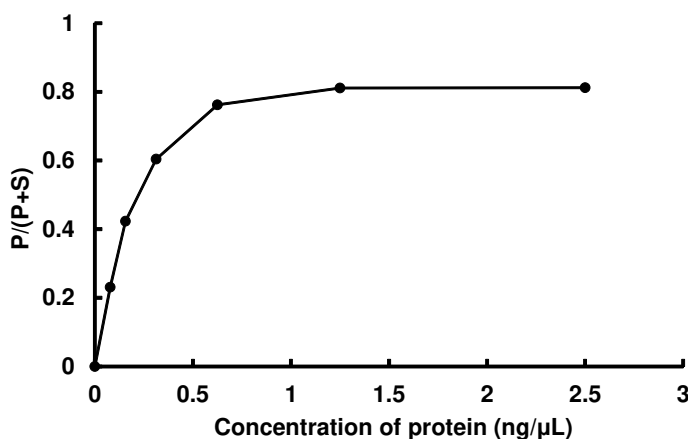
SDS-PAGE



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 μM