

## Product Information

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

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

#### Product description

Full-length human AMPK $\alpha$ 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-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.

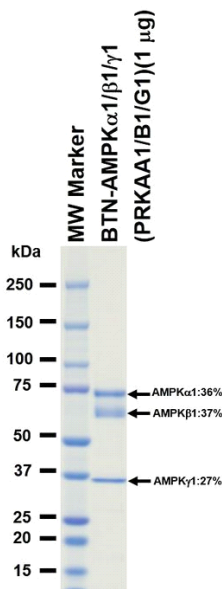
#### Storage buffer:

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.

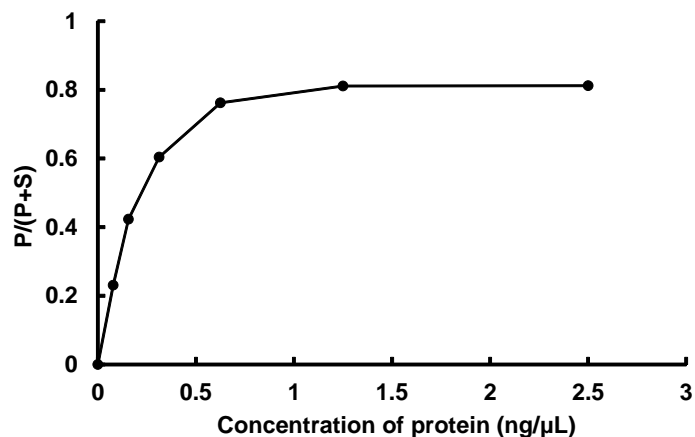
#### SDS-PAGE



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  $\mu$ M