

## Product Information

### AMPK $\alpha$ 2/ $\beta$ 2/ $\gamma$ 1(PRKAA2/B2/G1)

Product Number : 02-148

#### Product description

Full-length human AMPK $\alpha$ 2 [1-552(end) amino acids of accession number NP\_006243.2] was co-expressed as N-terminal GST-fusion protein (89 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. GST-AMPK $\alpha$ 2/ $\beta$ 2/ $\gamma$ 1 was purified by using glutathione sepharose chromatography and activated with His-tagged CaMKK1. Activated GST-AMPK $\alpha$ 2/ $\beta$ 2/ $\gamma$ 1 was purified by using glutathione sepharose chromatography.

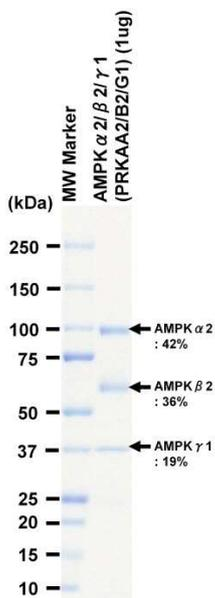
#### Storage buffer:

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.

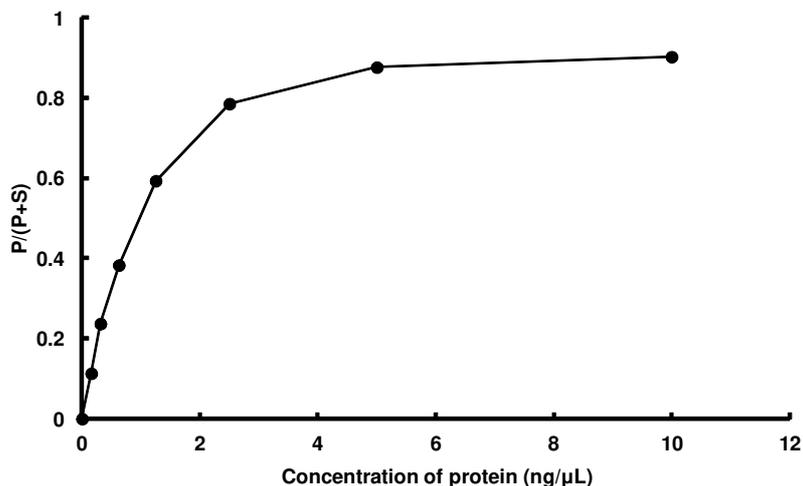
#### SDS-PAGE



Purity: 97 %

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