

## Data Sheet

### AMPK (A1/B1/G1), His-tag

Human, recombinant, C-terminal His-tag  
**Catalog #:** 40025  
**Lot #:** 190613-1      **Conc.:** 0.30 mg/ml

**Fomulated in:** 50 mM sodium phosphate, pH 7.0, 300 mM NaCl, 150 mM imidazole, 0.25 mM DTT, 25% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

**References:**

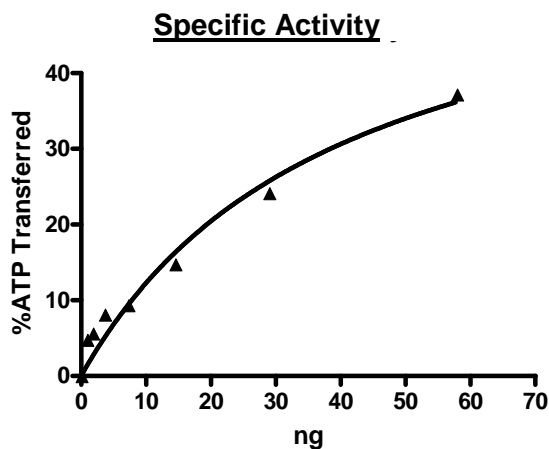
1. Minokoshi, Y. *et al. Nature* **428**: 569-574, 2004.
2. Hardie, D G. *et al. Eur J Biochem.* 1997 Jun 1; **246(2)**:259-73.

**Description:** Recombinant full-length human AMPK (combination of A1/B1/G1 subunits) was expressed by baculovirus in Sf9 insect cells using a C-terminal His-tag. The Genbank accession number for the three subunits (A1/B1/G1) is NM\_006251, NM\_006253, and NM\_002733, respectively. This protein has been activated with CAMKK.

**Specific Activity:**  $\geq 300$  pmol/min/ $\mu\text{g}$   
 Assay Buffer: Assay was done in Kinase buffer containing 1 mM DTT using AMARA peptide as a substrate (0.1 mg/ml), 20  $\mu\text{M}$  ATP and 100  $\mu\text{M}$  AMP. Reaction was done at  $30^{\circ}\text{C}$  for 45 min. Amount of ATP transferred was calculated using Kinase-Glo reagent (Promega)

**Application:** Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

### Quality Assurance



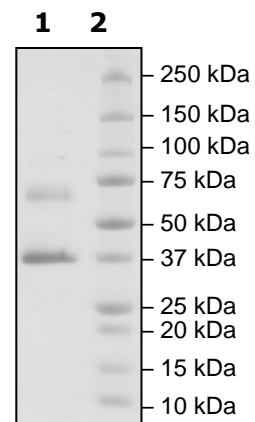
**4-20% SDS-PAGE  
 Coomassie staining**

**Lane 1:**  
AMPK

**Lane 2:**  
Protein Marker

**Purity:**  $\geq 90\%$

**MW:**  
 AMPKa1: 65kDa  
 AMPKg1: 38kDa  
 AMPKb1: 31kDa



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.259.5734**  
 Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)