

## Data Sheet

### AMPK (A2/B1/G1)

Human, recombinant, C-terminal His-tag

**Catalog #:** 40024

**Lot #:** 141008

**Conc.:** 0.1 mg/ml

**Formulated in:** 50 mM sodium phosphate, pH 7.0, 300 mM NaCl, 150 mM imidazole, 0.1 mM PMSF, 0.25 mM DTT, and 25% glycerol.

**Stability:** >6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

**References:**

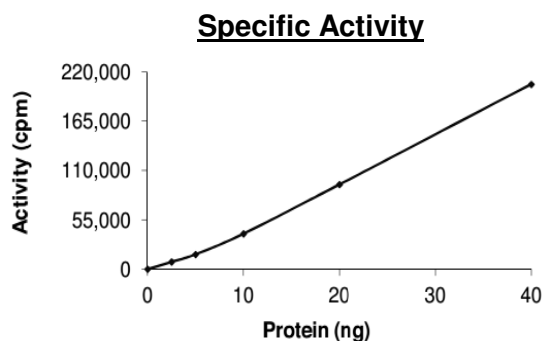
1. Viollet, B., *Biochem Soc Trans.* 2003;**31**:216-219
2. Foretz, M., *Diabetes.* 2005;**54(5)**:1331-1339.

**Description:** Recombinant full-length human AMPK (combination of A2/B1/G1 subunits) was expressed by baculovirus in Sf9 insect cells using a C-terminal His-tag. The gene accession number for the three subunits (A2/B1/G1) is NM\_006252, NM\_006253, and NM\_002733.

**Specific Activity:** 280 pmol/min/ $\mu\text{g}$   
 Assay Buffer: 5 mM MOPS, 2.5 mM  $\beta$ -glycerophosphate, 5 mM  $\text{MgCl}_2$ , 1 mM EGTA, 0.4 mM EDTA, 0.05 mM DTT, 2 mM ATP. Incubate AMPK A2/B1/G1 with 0.2 mg/ml SAMStide peptide substrate, 0.1 mM AMP, and  $[^{33}\text{P}]$ -ATP at  $30^{\circ}\text{C}$  for 15 minutes, then spot reaction on phosphocellulose paper, fix in 1% phosphoric acid, and assay with a scintillation counter.

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

### Quality Assurance



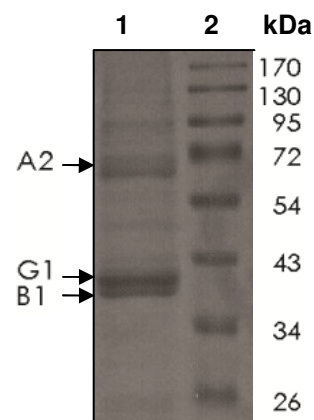
**10% SDS-PAGE  
 Coomassie staining**

**Lane 1:**  
AMPK A2 →

**Lane 2:**  
Protein Marker

**MW:**  
69 kDa (A2)  
38 kDa (B1)  
40 kDa (G1)

**Purity:**  $\geq 80\%$



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

To place your order, please contact us by Phone **1.858.481.8694** 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)