

## AMPK (A1/B2/G1), Active

Full-length recombinant protein expressed in Sf9 cells

#### Catalog # P50-10H Lot # Q224-1

## **Product Description**

Recombinant full-length human AMPK (combination of A1/B2/G1 subunits) was expressed by baculovirus in Sf9 insect cells using C-terminal His tags. The gene accession numbers for the three subunits (A1/B2/G1) are <u>NM 006251</u>, <u>NM 005399</u>, and <u>NM 002733</u>.

#### **Gene Aliases**

Subunit A1: PRKAA1, MGC33776, MGC57364 Subunit B2: PRKAB2, MGC61468 Subunit G1: PRKAG1, AMPKG, MGC8666

## **Formulation**

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.2mM DTT, 25% glycerol.

## Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

## **Scientific Background**

AMPK (A1/B2/G1) is a member of the AMPK family which are heterotrimeric proteins consisting of an alpha catalytic subunit, and non-catalytic beta and gamma subunits. AMPKs are an important energy-sensing enzyme group in the cells that monitor energy status particularly in response to stress (1). AMPKs regulate fatty acid and cholesterol synthesis by regulating the key rate-limiting enzymes acetyl-CoA carboxylase and hydroxy betamethylglutaryl-CoA reductase. The  $\beta$  subunit may be a positive regulator of AMPK activity and is highly expressed in skeletal muscle (2).

## References

- 1 Viollet, B. et al: Physiological role of AMP-activated protein kinase (AMPK): insights from knockout mouse models. Biochem. Soc. Trans. 2003; 31; 216-219.
- Thornton, C. et al: Identification of a novel AMP-activated 2. protein kinase beta subunit isoform that is highly expressed in skeletal muscle. J. Biol. Chem. 273: 12443-12450, 1998.

#### To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com www.signalchem.com

## **Specific Activity**



The specific activity of AMPK was determined to be 390 nmol /min/mg as per activity assay protocol.

## **Purity**



The purity of AMPK was determined to be >70% by densitometry, approx. MW ~68kDa (A1), ~36kDa (B2), and ~40kDa (G1).

## AMPK (A1/B2/G1), Active

Full-length recombinant protein expressed in Sf9 cells

Catalog Number Specific Activity Lot Specific Number Purity Concentration Stability Storage & Shipping

P50-10H 390 nmol/min/mg Q224-1

>70% 0.1 μg/μl 1yr At -70°C from date of shipment

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

# Activity Assay Protocol

**Reaction Components** 

#### Active Kinase (Catalog #: P50-10H)

Active AMPK ( $0.1\mu g/\mu l$ ) diluted with Kinase Dilution Buffer III (Catalog #: K23-09) and assayed as outlined in sample activity plot. (Note: these are suggested working dilutions and it is recommended that the researcher perform a serial dilution of Active AMPK for optimal results).

**Kinase Dilution Buffer III** (Catalog #: K23-09)

Kinase Assay Buffer I (Catalog #: K01-09) diluted at a 1:4 ratio (5X dilution) with 50ng/µl BSA.

Kinase Assay Buffer I (Catalog #: K01-09)

Buffer components: 25mM MOPS, pH 7. 2, 12.5mM  $\beta$ -glycerol-phosphate, 25mM MgCl<sub>2</sub>, 5mM EGTA, 2mM EDTA. Add 0.25mM DTT to Kinase Assay Buffer prior to use.

## [<sup>33</sup>P]-ATP Assay Cocktail

Prepare 250µM [<sup>33</sup>P]-ATP Assay Cocktail in a designated radioactive working area by adding the following components: 150µl of 10mM ATP Stock Solution (Catalog #: A50-09), 100µl [<sup>33P</sup>]-ATP (1mCi/100µl), 5.75ml of Kinase Assay Buffer I (Catalog #: K01-09). Store 1ml aliquots at -20°C.

10mM ATP Stock Solution (Catalog #: A50-09)

Prepare ATP stock solution by dissolving 55mg of ATP in 10ml of Kinase Assay Buffer I (Catalog #: K01-09). Store  $200\mu$ l aliquots at -20°C.

Substrate (Catalog #: S07-58)

SAMS synthetic peptide substrate (HMRSAMSGLHLVKRR) diluted in distilled H<sub>2</sub>O to a final concentration of 1mg/ml.

#### Assay Protocol

- Step 1. Thaw [33P]-ATP Assay Cocktail in shielded container in a designated radioactive working area.
- Step 2. Thaw the Active AMPK, Kinase Assay Buffer, Substrate and Kinase Dilution Buffer on ice.
- Step 3. In a pre-cooled microfuge tube, add the following reaction components bringing the initial reaction volume up to 20μl:
  - Component 1. 10µl of diluted Active AMPK (Catalog # P50-10H)
  - **Component 2.** 5µl of 1mg/ml stock solution of substrate (Catalog # \$07-58)

**Component 3.** 5µl of 0.5mM AMP solution (Catalog # A46-09)

- Step 4. Set up the blank control as outlined in step 3, excluding the addition of the substrate. Replace the substrate with an equal volume of distilled H<sub>2</sub>O.
- Step 5. Initiate the reaction by the addition of 5 μl [<sup>33</sup>P]-ATP Assay Cocktail bringing the final volume up to 25μl and incubate the mixture in a water bath at 30°C for 15 minutes.
- **Step 6.** After the 15 minute incubation period, terminate the reaction by spotting 20 μl of the reaction mixture onto individual pre-cut strips of phosphocellulose P81 paper.
- Step 7. Air dry the pre-cut P81 strip and sequentially wash in a 1% phosphoric acid solution (dilute 10ml of phosphoric acid and make a 1L solution with distilled H<sub>2</sub>O) with constant gentle stirring. It is recommended that the strips be washed a total of 3 intervals for approximately 10 minutes each.
- Step 8. Count the radioactivity on the P81 paper in the presence of scintillation fluid in a scintillation counter.
- Step 9. Determine the corrected cpm by removing the blank control value (see Step 4) for each sample and calculate the kinase specific activity as outlined below.

## Calculation of [P<sup>33</sup>]-ATP Specific Activity (SA) (cpm/pmol)

Specific activity (SA) = cpm for 5 µl [<sup>33</sup>P]-ATP / pmoles of ATP (in 5 µl of a 250 µM ATP stock solution, i.e., 1250 pmoles)

#### Kinase Specific Activity (SA) (pmol/min/µg or nmol/min/mg)

Corrected cpm from reaction / [(SA of <sup>33</sup>P-ATP in cpm/pmol)\*(Reaction time in min)\*(Enzyme amount in µg or mg)]\*[(Reaction Volume) / (Spot Volume)]

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: <u>orders@signalchem.com</u> <u>www.signalchem.com</u>

#### FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.