

AMPKG3 Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog # A60-30CG

Lot # T848-3

Product Description

Recombinant full-length human AMPKG3 subunit was expressed by baculovirus in Sf9 insect cells using the N-terminal GST and C-terminal His tags. The gene accession numbers is [NM_017431](#).

Gene Aliases

PRKAG3

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 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.

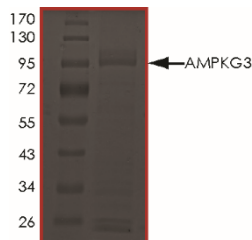
Scientific Background

5'-AMP-activated protein kinase subunit gamma-3 (AMPKG3) is an AMP/ATP-binding subunit of AMP-activated protein kinase (AMPK), an energy sensor protein kinase that plays a key role in regulating cellular energy metabolism. Ablation of the γ 3 subunit is associated with impaired glycogen resynthesis in skeletal muscle. In contrast, naturally occurring mutations in the γ 3 subunit increase glycogen storage in skeletal muscle, which may enhance glycogen storage by shifting the metabolic fate of glucose from oxidation to storage.

References

1. Barnes BR, et al. 5'-AMP-activated protein kinase regulates skeletal muscle glycogen content and ergogenics. *FASEB J.* 19:773-9, 2005.
2. Mu J, et al. A role for AMP-activated protein kinase in contraction- and hypoxia-regulated glucose transport in skeletal muscle. *Mol Cell.* 7:1085-94, 2001.

Purity



The purity of AMPKG3 was determined to be **>80%** by densitometry, approx. MW **96~108kDa**.

AMPKG3 Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog #	A60-30CG
Lot #	T848-3
Purity	>80%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	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.

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

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