# AMPK alpha 1 (T172D) [GST-tagged]

**Kinase** 

Alternate Names: 5'-AMP-activated protein kinase catalytic subunit alpha-1, Acetyl-CoA carboxylase kinase, Hydroxymethylglutaryl-CoA reductase kinase, HMGCR kinase, Tau-protein kinase PRKAA1

Cat. No.	66-0040-050
Lot. No.	30319

Quantity: 50 µg Storage: -70°C

FOR RESEARCH USE ONLY

NOT FOR USE IN HUMANS



# **CERTIFICATE OF ANALYSIS Page 1 of 2**

Protein Sequence: Please see page 2

### Background

Protein ubiquitylation and protein phosphorylation are the two major mechanisms that regulate the functions of proteins in eukaryotic cells. However, these different posttranslational modifications do not operate independently of one another, but are frequently interlinked to enable biological processes to be controlled in a more complex and sophisticated manner. Studying how protein phosphorylation events control the ubiquitin system and how ubiguitylation regulates protein phosphorylation has become a focal point of the study of cell regulation and human disease. Cloning of human 5'-AMP-activated protein kinase catalytic subunit alpha-1 (AMPK alpha 1) was first described by Stapleton et al. (1996). An example of such interplay between phosphorylation and ubiquitylation has been highlighted in recent studies indicating that AMPK alpha, along with AMPK kinases NUAK1 and MARK4, can be ubiquitylated with atypical ubiguitin chains. The deubiguitylating enzyme (DUB) found to remove these ubiquitin chains from both NUAK1 and MARK4 has been identified as USP9X. AMPK activation has also been shown to increase the expression of the E3 ubiquitin ligases MAFBx/Atrogin-1 and MuRF1. These ubiquitin ligases regulate key cardiac transcription factors to control cardiomyocyte mass and remodeling, thus suggesting another mechanism by which AMPK may function in the heart. The relevance of AMPK ubiquitylation in cardiac disease has yet to be



Species: rat

Source: E. coli

Quantity: 50 µg

Concentration: 2.0 mg/ml

Formulation: 50 mM Tris/HCl pH7.5, 0.1 mM EGTA, 150 mM NaCl, 0.1% ß-Mercaptoethanol, 270 mM sucrose, 0.03% Brij-35, 1 mM Benzamidine, 0.2 mM PMSF

Molecular Weight: ~64.6 kDa

Purity: >95% by InstantBlue™ SDS-PAGE

Stability/Storage: 12 months at -70°C; aliquot as required

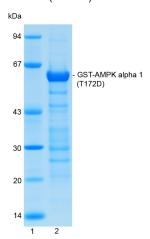
## **Quality Assurance**

### **Purity:**

**ORDERS / SALES SUPPORT** 

Email: sales.support@ubiquigent.com

4-12% gradient SDS-PAGE InstantBlue<sup>™</sup> staining Lane 1: MW markers Lane 2: 2.5 µg GST-AMPK alpha 1 (T172D)



### **Protein Identification:**

Confirmed by mass spectrometry.

### Activity Assay:

The specific activity of GST-AMPK alpha 1 (T172D) was determined using the method described by Hastie et al. (2006) with the enzyme being assayed at several concentrations. GST-AMPK alpha 1 (T172D) was incubated for 10 minutes at 30°C in kinase reaction buffer in the presence of AMARA peptide substrate (300 µM) and [y-32P]ATP (100 µM). Duplicate reactions were stopped by spotting the assay mixture onto Whatman P81 paper - capturing the phosphorylated substrate. The radioactivity incorporated was measured on a scintillation counter and the enzyme's mean specific activity was calculated.

GST-AMPK alpha 1 (T172D) specific activity: 41.73 Units/mg (64.27 Units/ml)

1 Unit = 1 nmole of phosphate incorporated into the substrate in 1 minute

### Substrate: AMARA peptide (AMARAASAAALARRR)

(9AM-5PM UTC)

© Ubiquigent 2014. Unless otherwise noted, Ubiquigent, Ubiquigent logo and all other trademarks are the property of Ubiquigent, Ltd. **UK HQ and TECHNICAL SUPPORT** International: +44 (0) 1382 381147 (9AM-5PM UTC)

Limited Terms of Use: For research use only. Not for use in humans or for diagnostics. Not for distribution or resale in any form, modification or derivative OR for use in providing services to a third party (e.g. screening or profiling) without the written permission of Ubiquigent, Ltd.

Email: tech.support@ubiquigent.com Email services@ubiquigent.com for enquiries regarding compound profiling and/or custom assay development services.

US/Canada: +1-617-245-0020

Lot-specific COA version tracker: v1.0.0

International: +1-617-245-0003 Ubiquigent US Toll-Free: 1-888-4E1E2E3 (1-888-431-3233) www.ubiguigent.com

Continued on page 2

Dundee, Scotland, UK

# AMPK alpha 1 (T172D) [GST-tagged]

Kinase

Alternate Names: 5'-AMP-activated protein kinase catalytic subunit alpha-1, Acetyl-CoA carboxylase kinase, Hydroxymethylglutaryl-CoA reductase kinase, HMGCR kinase, Tau-protein kinase PRKAA1



Quantity: 50 µg Storage: -70°C

NOT FOR USE IN HUMANS



**CERTIFICATE OF ANALYSIS Page 2 of 2** 

# Background

### Continued from page 1

tested directly, but it likely represents an important mechanism that occurs in common cardiac diseases that may be targeted for therapy (Zungu *et al.*, 2011).

### **References:**

Hastie CJ, McLauchlan HJ, Cohen P (2006) Assay of protein kinases using radiolabeled ATP: a protocol. *Nat Protoc* 1, 968-71.

Stapleton D, Mitchelhill KI, Gao G, Widmer J, Michell BJ, Teh T, et al. (1996) Mammalian AMP-activated protein kinase subfamily. J Biol Chem 271, 611-614.

Stein SC, Woods A, Jones NA, Davison MD, Carling D (2000) The regulation of AMP-activated protein kinase by phosphorylation. *Biochem J* **345**, 437-443.

Zungu M, Schisler JC, Essop MF, McCudden C, Patterson C and Willis MS (2011) Regulation of AMPK by the ubiquitin proteasome system. *Am J Pathol* **178**, 4-11.

**Physical Characteristics** 

### Continued from page 1

Protein Sequence:

**MSPILGYWKIKGLVQPTRLLLEYLEEKYEEH** LYERDEGDKWRNKKFELGLEFPNLPYYIDGD VKLTQSMAIIRYIADKHNMLGGCPKERAEISM LEGAVLDIRYGVSRIAYSKDFETLKVDFL SKLPEMLKMFEDRLCHKTYLNGDHVTHPD FMLYDALDVVLYMDPMCLDAFPKLVCFK **KRIEAIPQIDKYLKSSKYIAWPLQGWQATF** GGGDHPPKSDLVPRGSEFAMEQKLISEEDL GGG**E**KOKHDGRVKIGHYILGDTLGVGTFG **KVKVGKHELTGHKVAVKILNRQKIRSLDV** VGKIRREIONLKLFRHPHIIKLYOVISTPS DIFMVMEYVSGGELFDYICKNGRLDEKESR RLFQQILSGVDYCHRHMVVHRDLKPENVLL DAHMNAKIADFGLSNMMSDGEFLRDSCGSP NYAAPEVISGRLYAGPEVDIWSSGVILYALL CGTLPFDDDHVPTLFKKICDGIFYTPQYL NPSVISLLKHMLQVDPMKRATIKDIREHEW FKQDLPKYLFPEDPSYSSTMIDDEALKEVCEK FECSEEEVLRSITLAAARDRLD

Tag (**bold text**): N-terminal GST and N-terminal MYC Protease cleavage site: Thrombin (LVPR▼GS) AMPK alpha 1 (regular text): Start **bold italics** (amino acid residues 3-308). Non-AMPK sequence derived from the cloning vector (**bold**.

underlined). AMPK alpha 1 has a T172D mutation to mimic the activation of the enzyme through physiophysician of Thr172 by PDK1

of the enzyme through phosphorylation of Thr172 by PDK1 (Stein *et al.*, 2000). Accession number: AAC52355



### **ORDERS / SALES SUPPORT**

 International:
 +1-617-245-0003

 US Toll-Free:
 1-888-4E1E2E3 (1-888-431-3233)

 Email:
 sales.support@ubiquigent.com

#### **UK HQ and TECHNICAL SUPPORT**

 International:
 +44 (0) 1382 381147
 (9AM-5PM UTC)

 US/Canada:
 +1-617-245-0020
 (9AM-5PM UTC)

 Email:
 tech.support@ubiquigent.com

Email services@ubiquigent.com for enquiries regarding compound profiling and/or custom assay development services. © Ubiquigent 2014. Unless otherwise noted, Ubiquigent, Ubiquigent logo and all other trademarks are the property of Ubiquigent, Ltd.

Limited Terms of Use: For research use only. Not for use in humans or for diagnostics. Not for distribution or resale in any form, modification or derivative QR for use in providing services to a third party (e.g. screening or profiling) without the written permission of Ubiquigent, Ltd.

Lot-specific COA version tracker: v1.0.0