BRSK2 [6His-tagged]

Kinase

Alternate Names: BR Serine/Threonine Kinase 2; SAD1; PEN11B

Cat. No. 66-0001-050 Quantity: 50 μg **Lot. No. 2138** Storage: -70°C

FOR RESEARCH USE ONLY NOT FOR USE IN HUMANS



CERTIFICATE OF ANALYSIS Page 1 of 2

Background

by Sir Philip Cohen

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 ubiquitylation requlates protein phosphorylation has become a focal point of the study of cell regulation and human disease. Cloning of human Brain Specific Kinase 2 (BRSK2) was first described by Miura et al. (1998). BRSK2 is a member of the subfamily of protein kinases that include the AMP-activated protein kinase (AMPK) and, like AMPK itself, is activated by the tumour suppressor kinase LKB1 (Lizcano et al., 2004). As implied by its name BRSK2 is expressed in the brain where it plays an essential role in controlling neuronal cell polarisation. BRSK2 contains a ubiquitin-like domain adjacent to the kinase catalytic domain (Al-Hakim et al., 2008).

References:

Al-Hakim AK, Zagorska A, Chapman L, Deak M, Peggie M, Alessi DR (2008) Control of AMPK-related kinases by USP9X and atypical Lys(29)/Lys(33)-linked polyubiquitin chains. *Biochem J* 411, 249-60.

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

Continued on page 2

Physical Characteristics

Species: human **Protein Sequence:** Please see page 2

Source: Sf21 insect cell-baculovirus

expression

Quantity: 50 µg

Concentration: 1 mg/ml

Formulation: 50 mM Tris/HCl pH 7.5, 0.1 mM EGTA, 150 mM NaCl, 270 mM sucrose, 0.03% Brij, 0.1% β-Mercaptoethanol, 1 mM Benzamidine, 0.2 mM PMSF

Molecular Weight: ~78.3 kDa

Purity: >95% by InstantBlue™ SDS-PAGE

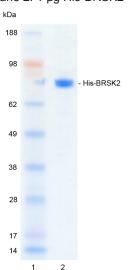
Stability/Storage: 12 months at -70°C;

aliquot as required

Quality Assurance

Purity:

4-12% gradient SDS-PAGE InstantBlue™ staining Lane 1: MW markers Lane 2: 1 µg His-BRSK2



Protein Identification:

Confirmed by mass spectrometry.

Activity Assay:

The specific activity of His-BRSK2 was determined using the method described by Hastie *et al.* (2006) with the enzyme being assayed at several concentrations. His-BRSK2 was incubated for 10 minutes at 30°C in kinase reaction buffer in the presence of CHKtide substrate (300µM) and [γ -3²P]ATP (100µM). 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.

His-BRSK2 specific activity:

333.7 Units/mg (333.7 Units/ml)

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

Substrate: CHKtide (KKKVSRSGLYRSPSMPENLNRPR)



Dundee, Scotland, UK

ORDERS / SALES SUPPORT

International: +1-617-245-0003 US Toll-Free: 1-888-4E1E2E3 (1-888-431-

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 2011**. 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 OR 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

BRSK2 [6His-tagged]

Kinase

Alternate Names: BR Serine/Threonine Kinase 2; SAD1; PEN11B

Cat. No. 66-0001-050 Quantity: 50 μg **Lot. No. 2138** Storage: -70°C

FOR RESEARCH USE ONLY NOT FOR USE IN HUMANS



CERTIFICATE OF ANALYSIS Page 2 of 2

Background

by Sir Philip Cohen

Continued from page 1

Miura K, Masuzaki H, Ishimaru T, Niikawa N, Jinno Y (1998) A Hhal/BstUI polymorphism in a novel gene at human chromosome 11p15.5. *J Hum Genet* **90**, 283-4.

Lizcano JM, Goransson O, et al. (2004) LKB1 is a master kinase that activates 13 kinases of the AMPK subfamily, including MARK/PAR-1. *EMBO J* 23, 833-43.

Background kindly written by:

Sir Philip Cohen FRS, FRSE University of Dundee

Director of the Medical Research Council Protein Phosphorylation Unit (1990-2012)

Director of the Scottish Institute for Cell Signalling incorporating the Protein Ubiquitylation Unit (2008-2012)

Co-Director of the Division of Signal Transduction Therapy (1998-2012)

Deputy Director of the Division of Signal Transduction Therapy (from July 2012)

Professor Cohen's research group is studying the interplay between protein phosphorylation and protein ubiquitylation in the regulation of innate immunity.

Physical Characteristics

Continued from page 1

Protein Sequence:

 ${\tt MSYYHHHHHHDYDIPTT} {\tt ENLYFQG} {\tt AMGS} {\tt \textit{T}}$ STGKDGGAQHAQYVGPYRLEKTLGKGQT GLVKLGVHCVTCQKVAIKIVNREKLS ESVLMKVEREIAILKLIEHPHVLKLHD VYENKKYLYLVLEHVSGGELFDYLVK KGRLTPKEARKFFRQIISALDFCHSH SICHRDLKPENLLLDEKNNIRIADFG MASLOVGDSLLETSCGSPHYACPEVIR GEKYDGRKADVWSCGVILFALLVGALPFD DDNLRQLLEKVKRGVFHMPHFIPPDC OSLLRGMIEVDAARRLTLEHIOKHIWYIG GKNEPEPEQPIPRKVQIRSLPSLEDIDP DVLDSMHSLGCFRDRNKLLQDLLSEEEN QEKMIYFLLLDRKERYPSQEDEDLPPR NEIDPPRKRVDSPMLNRHGKRRPERKS MEVLSVTDGGSPVPARRAIEMAQH GQRSRSISGASSGLSTSPLSSPRVT PHPSPRGSPLPTPKGTPVHTPKESPAGT PNPTPPSSPSVGGVPWRARLNSIKNS FLGSPRFHRRKLQVPTPEEMSNLTPESS PELAKKSWFGNFISLEKEEQIFVVIKDK PLSSIKADIVHAFLSIPSLSHSVISQTS FRAEYKATGGPAVFQKPVKFQVDI TYTEGGEAQKENGIYSVTFTLLSGPSR RFKRVVETIQAQLLSTHDPPAAQHLSEPP PPAPGLSWGAGLKGQKVATSYESSL

Tag (**bold text**): N-terminal His Protease cleavage site: TEV (<u>ENLYF ▼QG</u>) BRSK2 (regular text): Start **bold italics** (amino acid residues 2-674) Accession number: AAP97725.1



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 2011**. 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 OR 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