# C-Raf (Y340D; Y341D) [GST-tagged]

Kinase

Alternate Names: RAF proto-oncogene serine/threonine-protein kinase, Proto-oncogene c-RAF, RAF1

Cat. No. Lot. No.

66-0024-050 30303

Quantity: 50 µg Storage: -70°C

FOR RESEARCH USE ONLY

NOT FOR USE IN HUMANS



**CERTIFICATE OF ANALYSIS Page 1 of 2** 

## Background

There are three Raf kinase family members, all serine/threonine kinases, identified as: A-Raf, B-Raf and C-Raf (Rahman et al., 2013). C-Raf acts as a regulatory link between the membrane-associated Ras GTPases and the MAPK/ERK cascade, and this critical regulatory link functions as a switch determining cell fate decisions including proliferation, differentiation, apoptosis, survival and oncogenic transformation (Chen et al., 2001). Cloning of the C-Raf gene was first described by Rapp et al. (1983). Regulation is a highly complex process involving membrane recruitment, protein-protein interactions, dimerization, and phosphorylation/dephosphorylation events. Ras-GTP recruits C-Raf to the membrane, thereby promoting its activation. The inactive conformation of C-Raf is maintained by autoinhibitory interactions occurring between the N-terminal regulatory and the C-terminal catalytic domains and by the binding of a 14-3-3 protein that contacts two phosphorylation sites on the kinase (Abraham et al., 2000). Although C-Raf is capable of mutating into an oncogene in experimental settings, it is B-Raf that is the true major player in carcinogenesis in humans. Approximately 20% of all examined human tumour samples display a mutated B-Raf gene (Emuss et al., 2005). The overwhelming majority of these mutations involve the exchange of a single amino acid (V600E) that can mimic the activation loop phosphorylation and immediately render the ki-

#### Continued on page 2

Ubiquigent www.ubiguigent.com

## **Physical Characteristics**

Species: human

Source: baculovirus expression vector system

Quantity: 50 µg

Concentration: 0.48 mg/ml

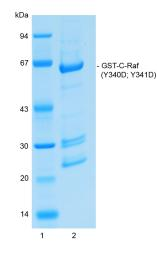
Formulation: 50 mM Tris/HCl pH7.5, 0.1 mM EGTA, 150 mM NaCl, 0.1% ß-Mercaptoethanol, 270 mM sucrose, 0.03% Brij-35, 1 mM Benzamidine, 0.2 mM PMSF

Molecular Weight: ~65.8 kDa

### **Quality Assurance**

#### **Purity:**

4-12% gradient SDS-PAGE InstantBlue<sup>™</sup> staining Lane 1: MW markers Lane 2: 2.5 µg GST-C-Raf (Y340D; Y341D)



Purity: >80% by InstantBlue™ SDS-PAGE

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

Protein Sequence: Please see page 2

# **Protein Identification:**

Confirmed by mass spectrometry.

#### Activity Assay:

The specific activity of GST-C-Raf (Y340D; Y341D) was determined using the method described by Hastie et al. (2006) with the enzyme being assayed at several concentrations. Initially, GST-C-Raf (Y340D; Y341D) (diluted in 50 mM Tris/HCl pH7.5, 0.1 mM EGTA, 1 mg/ml BSA, 10 mM DTT) was incubated with MKK1 (0.4 µg), p42MAPK (0.4 µg) and ATP (0.1 mM) in 50 mM Tris/ HCI pH7.5, 0.1 mM EGTA, 10 mM MgAc, 10 mM DTT buffer for 30 minutes at 30°C. A sample of this GST-C-Raf (Y340D; Y341D) reaction was then incubated for 10 minutes at 30°C in kinase reaction buffer in the presence of MBP substrate (0.33 mg/ml) 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-C-Raf (Y340D; Y341D) specific activity: 910028 Units/mg (436813 Units/ml)

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

Substrate: Myelin Basic Protein (MBP)

© Ubiquigent 2014. Unless otherwise noted, Ubiquigent, Ubiquigent logo and all other trademarks are the property of Ubiquigent, Ltd. **ORDERS / SALES SUPPORT UK HQ and TECHNICAL SUPPORT** International: +1-617-245-0003 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. US Toll-Free: 1-888-4E1E2E3 (1-888-431-3233) US/Canada: +1-617-245-0020 (9AM-5PM UTC) Email: sales.support@ubiquigent.com Email: tech.support@ubiquigent.com Email services@ubiquigent.com for enquiries regarding compound profiling and/or custom assay development services. Lot-specific COA version tracker: v1.0.0 Dundee, Scotland, UK

# C-Raf (Y340D; Y341D) [GST-tagged]

Kinase

Alternate Names: RAF proto-oncogene serine/threonine-protein kinase, Proto-oncogene c-RAF, RAF1

Cat. No. 66-0024-050 Lot. No. 30303

Quantity: 50 µg -70°C

Storage:

FOR RESEARCH USE ONLY

NOT FOR USE IN HUMANS



**CERTIFICATE OF ANALYSIS Page 2 of 2** 

### Background

# **Physical Characteristics**

Continued from page 1

#### Continued from page 1

nase domain fully active (Tran et al., 2005). Since B-Raf can also activate itself by homodimerisation and C-Raf by heterodimerisation, this mutation has a catastrophic effect by turning the ERK1/2 pathway constitutively active, driving an uncontrolled process of cell division (Garnett et al., 2005). Genetic mutations Y340D and Y341D have been shown to achieve maximum activation of C-Raf, simulating tyrosine phosphorylation and recruitment to the plasma membrane (Roy et al., 1997).

#### **References:**

Abraham D. Podar K. Pacher M. Kubicek M. Welzel N. Hemmings BA, et al. (2000) Raf-1-associated protein phosphatase 2A as a positive regulator of kinase activation. J Biol Chem 275, 22300-22304

Chen J, Fujii K, Zhang L, Roberts T and Fu H (2001) Raf-1 promotes cell survival by antagonizing apoptosis signal-regulating kinase 1 through a MEK-ERK independent mechanism. Proc Natl Acad Sci U S A 98, 7783-7788.

Emuss V. Garnett M. Mason C and Marais R (2005) Mutations of C-RAF are rare in human cancer because C-RAF has a low basa kinase activity compared with B-RAF. Cancer Res 65, 9719-9726.

Garnett MJ, Rana S, Paterson H, Barford D and Marais R (2005) Wild-type and mutant B-RAF activate C-RAF through distinct mechanisms involving heterodimerization. Mol Cell 20, 963-969.

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

Rahman MA, Salajegheh A, Smith RA and Lam AK (2013) B-Raf mutation: a key player in molecular biology of cancer. Exp Mol Pathol 95, 336-342.

Rapp UR, Goldsborough MD, Mark GE, Bonner TI, Groffen J, Reynolds FH, Jr., et al. (1983) Structure and biological activity of v raf, a unique oncogene transduced by a retrovirus. Proc Natl Acad Sci U S A 80, 4218-4222.

Roy S. Lane A. Yan J. McPherson R and Hancock JF (1997) Activity of plasma membrane-recruited Raf-1 is regulated by Ras via the Raf zinc finger. J Biol Chem 272, 20139-20145.

Tran NH, Wu X and Frost JA (2005) B-Raf and Raf-1 are regulated by distinct autoregulatory mechanisms. J Biol Chem 280, 16244-16253

#### **Protein Sequence: MSPILGYWKIKGLVQPTRLLLEYLEEKYEEH** LYERDEGDKWRNKKFELGLEFPNLPYYIDGD VKLTQSMAIIRYIADKHNMLGGCPKERAEISM LEGAVLDIRYGVSRIAYSKDFETLKVDFL SKLPEMLKMFEDRLCHKTYLNGDHVTHPD FMLYDALDVVLYMDPMCLDAFPKLVCFK **KRIEAIPQIDKYLKSSKYIAWPLQGWQAT** FGGGDHPPKSDLEVLFOGPLGSSOPKTPV PAORERAPVSGTOEKNKIRPRGORDSSD DWEIEASEVMLSTRIGSGSFGTVYKGKWHGD VAVKILKVVDPTPEOFOAFRNEVAVLRK TRHVNILLFMGYMTKDNLAIVTQWCEGSS LYKHLHVQETKFQMFQLIDIARQTAQGM DYLHAKNIIHRDMKSNNIFLHEGLTVKIGD FGLATVKSRWSGSQQVEQPTGSVLWMAPE VIRMODNNPFSFOSDVYSYGIVLYELMT GELPYSHINNRDQIIFMVGRGYASPDLSK LYKNCPKAMKRLVADCVKKVKEERPLFPQ ILSSIELLQHSLPKINRSASEPSLHRAAHTED INACTLTTSPRLPVF

Tag (bold text): N-terminal GST

Protease cleavage site: PreScission™ (LEVLFQ▼GP) C-Raf (regular text): Start bold italics (amino acid residues 306-648).

The enzyme has two mutations (Y340D and Y341D) to mimic phosphorylation and activation of C-Raf Accession number: NP 002871

Ubiquigent www.ubiquigent.com Dundee, Scotland, UK

**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 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