# **IKK epsilon** [GST-tagged]

Kinase

Alternate Names:	Inhibitor of nuclear factor kappa-B kinase subunit epsilon, I-kappa-B kinase epsilon, IKK-E, Inducible I kappa-B kinase, IKBKE, IKKI, KIAA0151

Cat. No.	66-0038-050
Lot. No.	30317

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. Inhibitor of IkB kinases (IKK) are key regulators of NF-KB signalling. Three IKK isoforms- $\alpha$ ,  $\beta$ , and ε-have been linked to oncogenesis (Hsu et al., 2012). IKK epsilon (IKKε) is a key regulator of innate immunity and a breast cancer oncogene, amplified in ~30% of breast cancers, that promotes malignant transformation through NF-κB activation (Zhou et al., 2013). Cloning of the IKK epsilon gene was first described by Shimada et al. (1999). IKK epsilon can be modified and regulated by K63-linked polyubiquitylation at lysine 30 and lysine 401. Tumour necrosis factor alpha (TNF $\alpha$ ) and interleukin-1ß (IL-1ß) stimulation can induce IKK epsilon K63-linked polyubiquitylation, and this modification is essential for IKK epsilon kinase activity, IKK epsilon-mediated NF-kB activation, and IKK epsilon-induced malignant transformation. Disruption of K63-linked ubiguitylation of IKK epsilon does not affect its overall struc-

### Continued on page 2



**ORDERS / SALES SUPPORT** International: +1-617-245-0003 US Toll-Free: 1-888-4E1E2E3 (1-888-431-3233) Email: sales.support@ubiquigent.com

Species: human

Source: baculovirus expression vector system

**Physical Characteristics** 

Quantity: 50 µg

Concentration: 0.34 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: ~107.3 kDa

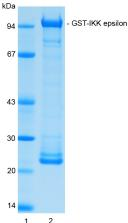
Purity: >60% by InstantBlue™ SDS-PAGE

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

## **Quality Assurance**

## **Purity:**

4-12% gradient SDS-PAGE InstantBlue<sup>™</sup> staining Lane 1: MW markers Lane 2: 2.5 µg GST-IKK epsilon



2

**Protein Identification:** 

Confirmed by mass spectrometry.

## Activity Assay:

The specific activity of GST-IKK epsilon was determined using the method described by Hastie et al. (2006) with the enzyme being assayed at several concentrations. GST-IKK epsilon was incubated for 10 minutes at 30°C in kinase reaction buffer in the presence of MBP substrate (1 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-IKK epsilon specific activity: 561.7 Units/mg (190.9 Units/ml)

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

Substrate: Myelin Basic Protein (MBP)

**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

# **IKK epsilon** [GST-tagged]

Kinase

Alternate Names:	Inhibitor of nuclear factor kappa-B kinase subunit epsilon, I-kappa-B kinase epsilon, IKK-E,
	Inducible I kappa-B kinase, IKBKE, IKKI, KIAA0151

Cat. No.	66-0038-050
Cat. No. Lot. No.	30317
FOR RESEARCH USE ONLY	

Quantity: 50 µg Storage: -70°C

NOT FOR USE IN HUMANS



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

## Background

### Continued from page 1

ture but impairs the recruitment of canonical NF- $\kappa$ B proteins. The ubiquitin E3 ligase complex involved in binding to and ubiquitylating IKK epsilon is cIAP1/cIAP2/TRAF2 (Zhou *et al.*, 2013).

### **References:**

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

Hsu S, Kim M, Hernandez L, Grajales V, Noonan A, Anver M, et al. (2012) IKK-epsilon coordinates invasion and metastasis of ovarian cancer. *Cancer Res* **72**, 5494-5504.

Shimada T, Kawai T, Takeda K, Matsumoto M, Inoue J, Tatsumi Y, et al. (1999) IKK-i, a novel lipopolysaccharide-inducible kinase that is related to IkappaB kinases. Int Immunol **11**, 1357-1362.

Zhou AY, Shen RR, Kim E, Lock YJ, Xu M, Chen ZJ, *et al.* (2013) IKKepsilon-mediated tumorigenesis requires K63-linked polyubiquitination by a cIAP1/cIAP2/TRAF2 E3 ubiquitin ligase complex. *Cell Rep* **3**, 724-733.

## **Physical Characteristics**

#### Continued from page 1

**Protein Sequence: MSPILGYWKIKGLVQPTRLLLEYLEEKYEEH** LYERDEGDKWRNKKFELGLEFPNLPYY IDGDVKLTQSMAIIRYIADKHNMLGGCP **KERAEISMLEGAVLDIRYGVSRIAYSKD** FETLKVDFLSKLPEMLKMFEDRLCHKTYLNGD HVTHPDFMLYDALDVVLYMDPMCLDAFP **KLVCFKKRIEAIPQIDKYLKSSKYIAWPLQG** WOATFGGGDHPPKSDLEVLFOGPLGSMOSTA NYLWHTDDLLGOGATASVYKARNKKSGEL VAVKVFNTTSYLRPREVQVREFEVLRKLN HONIVKLFAVEETGGSROKVLVMEYCSSG SLLSVLESPENAFGLPEDEFLVVLRCVVAG MNHLRENGIVHRDIKPGNIMRLVGEEGQSI YKLTDFGAARELDDDEKFVSVYGTEEYLH PDMYERAVLRKPQQKAFGVTVDLWSIGVT LYHAATGSLPFIPFGGPRRNKEIMYRIT TEKPAGAIAGAQRRENGPLEWSYTLPITC QLSLGLQSQLVPILANILEVEQAKCWGFDQF FAETSDILQRVVVHVFSLSQAVLHHIYIHAHN TIAIFQEAVHKQTSVAPRHQEYLFEGHLCV LEPSVSAQHIAHTTASSPLTLFSTAIPKGLA FRDPALDVPKFVPKVDLQADYNTAKGVL GAGYQALRLARALLDGQELMFRGLHWVMEV LQATCRRTLEVARTSLLYLSSSLGTERFSS VAGTPEIQELKAAAELRSRLRTLAEVLSRC SQNITETQESLSSLNRELVKSRDQVHEDR SIQQIQCCLDKMNFIYKQFKKSRMRPGLGYNE EQIHKLDKVNFSHLAKRLLQVFQEECVQKY QASLVTHGKRMRVVHETRNHLRLVGCS VAACNTEAQGVQESLSKLLEELSHQLLQDRAK GAOASPPPIAPYPSPTRKDLLLHMOELCEGM KLLASDLLDNNRIIERLNRVPAPPDV

Tag (**bold text**): N-terminal GST Protease cleavage site: PreScission™ (LEVLFQ▼GP) IKK epsilon (regular text): Start **bold italics** (amino acid residues 1-716). Accession number: NP\_054721



## **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