

# SGK3 [GST-tagged]

## Kinase and Substrate

**Alternate Names:** Serine/threonine-protein kinase Sgk3, Cytokine-independent survival kinase, Serum/glucocorticoid-regulated kinase 3, CISK, SGKL

**Cat. No.** 66-0021-050

**Lot. No.** 30300

**Quantity:** 50 µg

**Storage:** -70°C

FOR RESEARCH USE ONLY

NOT FOR USE IN HUMANS



CERTIFICATE OF ANALYSIS Page 1 of 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 ubiquitylation regulates protein phosphorylation has become a focal point of the study of cell regulation and human disease. The serum- and glucocorticoid-inducible protein kinase (SGK) family is made up of three isoforms, SGK1, 2, and 3, that are phosphatidylinositol-3-kinase (PI3-K)-dependent, serine/threonine kinases, with similar substrate specificity to protein kinase B (PKB). Consequently, the SGK family also regulates similar cell processes to the PKB kinases, including cell proliferation and survival (Bruhn *et al.*, 2013). SGK3 functions in parallel to PKB downstream of PI3-K. It shares ~55% identity with PKB in the kinase domain and is also a direct substrate of 3-phosphoinositide-dependent kinase 1 (PDK1). In contrast to PKB, SGK3 lacks the pleckstrin homology domain but contains a phox (PX) homology domain that binds phosphatidylinositol 3'-monophosphate and targets SGK3 to the early endosome (Wang *et al.*, 2014). SGK3 is unique within the SGK family by containing an N-terminal PX

## Physical Characteristics

**Species:** human

**Source:** *E. coli*

**Quantity:** 50 µg

**Concentration:** 4.2 mg/ml

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

**Molecular Weight:** ~42.4 kDa

**Purity:** >95% by InstantBlue™ SDS-PAGE

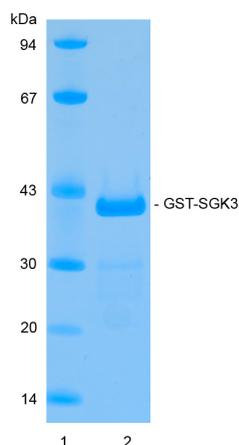
**Stability/Storage:** 12 months at -70°C; aliquot as required

**Protein Sequence:** Please see page 2

## Quality Assurance

**Purity:**  
4-12% gradient SDS-PAGE  
InstantBlue™ staining  
Lane 1: MW markers  
Lane 2: 2.5 µg GST-SGK3

**Protein Identification:**  
Confirmed by mass spectrometry.



Continued on page 2



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

# SGK3 [GST-tagged]

## Kinase and Substrate

**Alternate Names:** Serine/threonine-protein kinase Sgk3, Cytokine-independent survival kinase, Serum/glucocorticoid-regulated kinase 3, CISK, SGKL

**Cat. No.** 66-0021-050

**Lot. No.** 30300

**Quantity:** 50 µg

**Storage:** -70°C

FOR RESEARCH USE ONLY

NOT FOR USE IN HUMANS



CERTIFICATE OF ANALYSIS Page 2 of 2

### Background

Continued from page 1

domain which is shown to be important for targeting SGK3 to vesicle-like structures (Bruhn *et al.*, 2013). Cloning of the SGK3 gene was first described by Kobayashi *et al.* (1999). SGK3 (or SGK1) overexpression has been shown to increase phosphorylation of the ubiquitin E3 ligase Nedd4-2, which is known to inhibit Nedd4-2 activity (Lamothe and Zhang, 2013). Activated SGK3 can also promote estrogen/estrogen receptor (ER) dependent transcription and cell survival. Recent studies have established a clinical link between SGK3 and ER, underlining the importance of incorporating SGK3 as a new component in the assessment of breast cancer (Xu *et al.*, 2012).

### References:

Bruhn MA, Pearson RB, Hannan RD and Sheppard KE (2013) AKT-independent PI3-K signaling in cancer - emerging role for SGK3. *Cancer Manag Res* 5, 281-292.

Kobayashi T, Deak M, Morrice N and Cohen P (1999) Characterization of the structure and regulation of two novel isoforms of serum- and glucocorticoid-induced protein kinase. *The Biochemical Journal* 344 Pt 1, 189-197.

Lamothe SM and Zhang S (2013) The serum- and glucocorticoid-inducible kinases SGK1 and SGK3 regulate hERG channel expression via ubiquitin ligase Nedd4-2 and GTPase Rab11. *J Biol Chem* 288, 15075-15084.

Wang Y, Xu W, Zhou D, Neckers L and Chen S (2014) Coordinated regulation of serum- and glucocorticoid-inducible kinase 3 by a C-terminal hydrophobic motif and Hsp90-Cdc37 chaperone complex. *J Biol Chem* 289, 4815-4826.

Xu J, Wan M, He Q, Bassett RL, Jr., Fu X, Chen AC, *et al.* (2012) SGK3 is associated with estrogen receptor expression in breast cancer. *Breast Cancer Res Treat* 134, 531-541.

### Physical Characteristics

Continued from page 1

#### Protein Sequence:

**MSPILGYWKIKGLVQPTRLLLEYLEEKYEEH**  
**LYERDEGDKWRNKKFELGLEFPNLPYYIDGD**  
**VKLTQSMAIRYIADKHNMLGGCPKERAEISM**  
**LEGAVLDIRYGVSR IAYSKDFETLKVDVFL**  
**SKLPEMLKMFEDRLCHKTYLNGDHVTHPD**  
**FMLYDALDVVLYMDPMCLDAFPKLVCFK**  
**KRIEAIPOIDKYLKSSKYIAWPLQGWQAT**  
**FGGGDHPPKSD**LEVLFQGPLGSMQRDHTM  
DYKESCPSVSI PSSDEHREKKRFTVYKVLVS  
VGRSEWFVFRRYAEFDKLYNTLKKQFPAMAL  
KIPAKRIFGDNFDPDFIKQRRAGLNEFIQN  
LVRYPELYNHPDVRAFLQMDSPKHQSDPSE

Tag (**bold text**): N-terminal GST

Protease cleavage site: PreScission™ (LEVLFQ▼GP)

SGK3 (regular text): Start **bold italics** (amino acid residues 1-130).

Accession number: NP\_037389



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