

Certificate of Analysis



SGK3

Serum/glucocorticoid regulated kinase-like

Recombinant Human Active Protein Kinase

HGNC Symbol: SGK3

Synonyms: CISK, SGKL, SGK2

Product No.: 0198-0000-2

Lot: 004

Description: Human SGK3, full length, amino acids M₁-L₄₉₆ (as in NCBI/Protein entry NP_037389.4), activated, N-terminal GST-fusion protein, expressed in Sf9 insect cells

Product identity: SGK3 Lot 004, was confirmed as SGK3 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 83,186 Da

Expression: Baculovirus infected Sf9 cells

Purification: GST-Affinity Chromatography

Activation: With PDK1 (GenBank entry NM_002613)

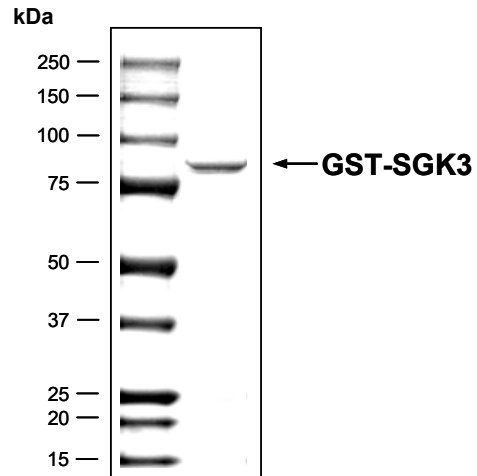
Storage buffer: 50 mM Tris-HCl pH 8.0, 100 mM NaCl, 5 mM DTT, 4 mM reduced glutathione, 20% glycerol

Storage temperature: -80°C
Avoid repeated freeze-thaw cycles!

Protein concentration: 0.201 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

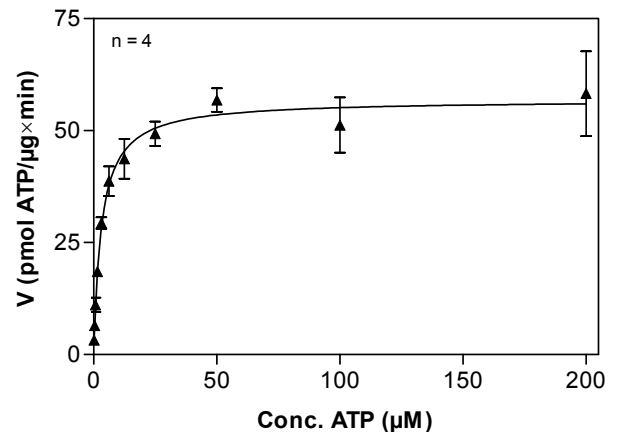
Biochemical Parameters:
Specific kinase activity (P_i transfer): 57 pmol/µg×min
ATP-K_M: 3.1 µM

**SGK3 Lot 004:
Coomassie stain**



2.0 µg GST-SGK3


**SGK3 Lot 004:
Determination of V_{max} and K_M value for ATP**



Determination of K_M value & Specific activity:

- Assay conditions:
 - 60 mM HEPES-NaOH, pH 7.5
 - 3 mM MgCl₂
 - 3 mM MnCl₂
 - 3 µM Na-orthovanadate
 - 1.2 mM DTT
 - 50 µg / ml PEG_{20,000}
 - ATP (variable)
 - Substrate: GSK3-derived peptide (R11-SGRARTSSFAEPGGK), 100 µg / ml
 - SGK3: 4.0 µg / ml
- Filter binding assay
MSPH membrane (Millipore)

Additional assay technology: SGK3 Lot 004

was also successfully tested by ProQinase for the use with the ADP-Glo™ Kinase assay from 

This product was manufactured at ProQinase in Freiburg, Germany and is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.

CoA V.1.1 SGK3_Lot004_V4.doc

SGK3

Product No.: 0198-0000-2

SGK3 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSR	YSKDFETLKV	120
121	DFLSKLP	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI	PQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PGIQ MQRDHT	240
241	SIPSSDEHRE	KKKRFTVYKV	LVSVGRSEWF	VFRRYA	EFDK	LYNTLKKQFP	300
301	IFGDNFDPDF	IKQRRAGLNE	FIQNLVRYPE	LYNHPDVRAF	LQMDSPKHQS	D	360
361	SQKLHSTSON	INLGPSGNPH	AKPTDFDFLK	VIGKGSFGKV	LLAKRKL	DGK	420
421	IVLNRKEQKH	IMAERNVLLK	NVKHPFLVGL	HYSFQTTEKL	YFVLDFVNGG	ELFFHLQ	480
481	SFPEHRARFY	AAEIASALGY	LHSIKIVYRD	LKPENILLDS	VGHVVL	TDFG	540
541	TTTTFCGTPE	YLAP	EVIKQ	YDNTVDW	WLGAVLYEMLY	GLPPFYCRDV	600
601	PLSLRPGVSL	TAWSILEELL	EKDRQNLGA	KEDFLEIQNH	PFESLSWAD	LVQKKIP	660
661	NPVAGPDDI	RNFDTAFTEE	TVPYSVCVSS	DYSIVNASVL	EADDAFVGFS	YAPPS	720

1-218: GST blue: SGK3 boxed: variation from RefSeq

SGK3 wt ¹ amino acid sequence							
1	MQRDHTMDYK	ESCPSVSIPS	SDEHREKKKR	FTVYKVLVSV	GRSEWVFVRR	YAEVDKLYNT	60
61	LKKQFPAMAL	KIPAKRIFGD	NFPDFIKQR	RAGLNEFIQN	LVRYPELYNH	PDVRAFLQMD	120
121	SPKHQSGPSE	DEDERSSQKL	HSTSONINLG	PSGNPHAKPT	DFDFLKVIGK	GSFGKVLAK	180
181	RKLDGK	VYAV	KVLQKKIVLN	RKEQKHIMAE	RNVLLKNVKH	PFLVGLHYSF	240
241	DFVNGGELFF	HLQERSFPE	HRARFYAAEI	ASALGYLHSI	KIVYRDLKPE	NILVDSVGHV	300
301	VLTDGFLCKE	GIAISDTTTT	FCGTPEYLAP	EVIRKQPYDN	TVDWWCLGAV	LYEMLYGLPP	360
361	FYCRDVAEMY	DNILHKPLSL	RPGVSLRAWS	ILEELLEKDR	QNLGAKEDF	LEIQNHPPFE	420
421	SLSWADLVQK	KIPPPFNPNV	AGPDDIRNFD	TAFTEETVPY	SVCVSSDYSI	VNASVLEADD	480
481	AFVGFSYAPP	SEDLFL					540

blue: SGK3 sequence expressed in fusionprotein Red: variant in fusionprotein

¹NCBI/Protein accession number NP_037389.4
G127D; V187F: documented sequence conflict, see UniProt Q96BR1/ENA entry AAF27051

This product was manufactured at ProQinase in Freiburg, Germany and is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.