

ProQinase™ SGK3

serum/glucocorticoid regulated kinase family member 3

Recombinant Human Active Protein Kinase

HGNC Symbol: SGK3

Synonyms: CISK, SGK1, 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 host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: With PDK1

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

For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. 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:

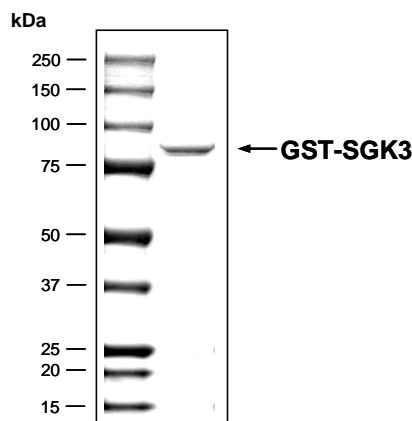
Specific kinase activity (P_i transfer): 57 pmol/µg × min

ATP-K_M: 3.1 µM

Additional assay technology:

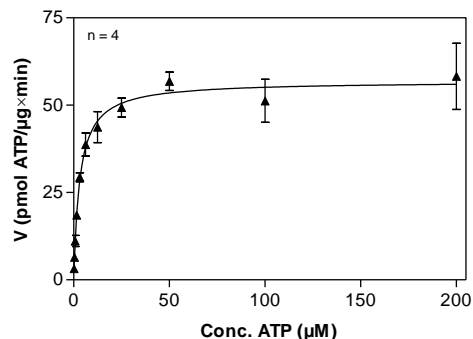
SGK3 Lot 004 was also successfully tested by Reaction Biology for the use with the ADP-Glo™ Kinase assay from Promega ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

**SGK3 Lot 004:
Coomassie stain**



2.0 µg GST-SGK3

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



- 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, 100 µg/ml
Kinase: 4 µg/ml
- Filter binding assay
MSPH membrane (Millipore)

ProQinase™ SGK3

Product No.: 0198-0000-2

GST-SGK3 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLIERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWAQTF	GGGDHPPKSD	PGIQMQRDHT	MDYKESCPSV	240
241	SIPSSDEHRE	KKKRFTVYKV	LVSVGRSEWF	VFRRYAEFDK	LYNTLKKQFP	AMALKIPAKR	300
301	IFGDNFDPDF	IKQRRAGLNE	FIQNLVRYPE	LYNHPDVRAF	LQMDSPKHQS	DPSEDEDEERS	360
361	SQKLHSTSQN	INLGPSGNPH	AKPTDFDFLK	VIGKGSFGKV	LLAKRKLDGK	FYAVKVLQKK	420
421	IVLNRKEQKH	IMAERNVLLK	NVKHPFLVGL	HYSFQTTEKL	YFVLDFVNGG	ELFFHLQREER	480
481	SFPEHRARFY	AAEIASALGY	LHSIKIVYRD	LKPENILLDS	VGHVLTDFG	LCKEGIAISD	540
541	TTTTFCGTPE	YLAPEVIRKQ	PYDNTVDWWC	LGAVLYEMLY	GLPPFYCRDV	AEMYDNILHK	600
601	PLSLRPGVSL	TAWSILEELL	EKDRQNLGA	KEDFLEIQNH	PFESLSWAD	LVQKKIPPPF	660
661	NPNVAGPDDI	RNFDTAFTEE	TVPYSVCVSS	DYSIVNASVL	EADDAFVGFS	YAPPSEDLFL	720

1-218: GST blue: SGK3

SGK3 wt ¹ Amino Acid Sequence							
1	MQRDHTMDYK	ESCPSVSIPS	SDEHREKKKR	FTVYKVLVSV	GRSEWFVFR	YAEFDKLYNT	60
61	LKKQFPAMAL	KIPAKRIFGD	NFDPDFIKQR	RAGLNEFIQN	LVRYPELYNH	PDVRAFLQMD	120
121	SPKHQSDPSE	DEDESSQKL	HSTSQNINLG	PSGNPHAKPT	DFDFLKVIGK	GSFGKVLLAK	180
181	RKLDGKFYAV	KVLQKKIVLN	RKEQKHIMAE	RNVLLKNVKH	PFLVGLHYSF	QTTEKLYFVL	240
241	DFVNGGELFF	HLQRERSFPE	HRARFYAAEI	ASALGYLHSI	KIVYRDLKPE	NILLDSVGHV	300
301	VLTFDGLCKE	GIAISDTTFT	FCGTPEYLAP	EVIRKQPYDN	TVDWWCLGAV	LYEMLYGLPP	360
361	FYCRDVAEMY	DNILHKPLSL	RPGVSLTAW	ILEELLEKDR	QNRLGAKEDF	LEIQNHPPFE	420
421	SLSWADLVQK	KIPPPFNPNV	AGPDDIRNFD	TAFTEETVPY	SVCVSSDYSI	VNASVLEADD	480
481	AFVGFYSYAPP	SEDLFL					540

blue: SGK3 sequence expressed in recombinant protein

¹[NCBI/Protein](https://www.ncbi.nlm.nih.gov/Protein/NP_037389.4) accession number NP_037389.4