

## ProQinase™ SGK1

serum/glucocorticoid regulated kinase 1

Recombinant Human Active Protein Kinase

HGNC Symbol: SGK1

Synonyms: SGK

Product No.: 0199-0000-2

Lot: 005

**Description:** Human SGK1, full length, amino acids M<sub>1</sub>-L<sub>431</sub> (as in [NCBI/Protein](#) entry NP\_005618.2), activated, N-terminal GST fusion protein, expressed in Sf9 insect cells

**Product identity:** SGK1 Lot 005 was confirmed as SGK1 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW**<sub>Fusion Protein</sub>: 74,950 Da

**Expression host:** Sf9 insect cells

**Purification:** GST-Affinity Chromatography

**Activation:** With PDK1

**Storage buffer:** 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 15 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:** 0.126 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

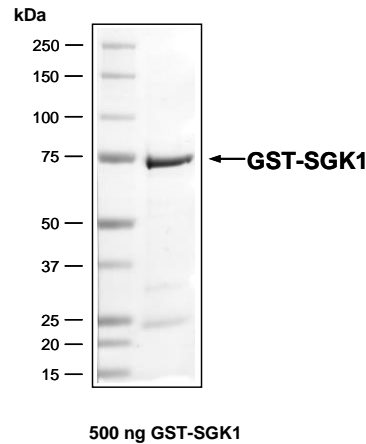
### Biochemical Parameters:

Specific kinase activity (P<sub>i</sub> transfer): 54 pmol/µg × min  
ATP-K<sub>M</sub>: 6.7 µM

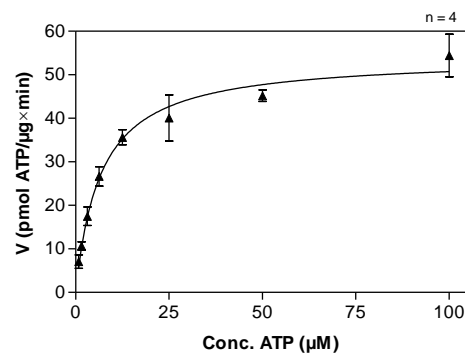
### Additional assay technology:

SGK1 Lot 005 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

### SGK1 Lot 005: Coomassie stain



### SGK1 Lot 005: Determination of V<sub>max</sub> and K<sub>M</sub> value for ATP



- Assay conditions:  
60 mM HEPES-NaOH, pH 7.5  
3 mM MgCl<sub>2</sub>  
3 mM MnCl<sub>2</sub>  
3 µM Na-orthovanadate  
1.2 mM DTT  
50 µg/ml PEG<sub>20,000</sub>  
ATP (variable)  
Substrate: GSK3-derived peptide, 40 µg/ml  
Kinase: 2 µg/ml
- Filter binding assay  
MSPH membrane (Millipore)

# ProQinase™ SGK1

Product No.: 0199-0000-2

GST-SGK1 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PGIQMTVKTE	<b>AAKGTLTYSR</b>	240
241	<b>MRGMVAILIA</b>	<b>FMKQRRMGLN</b>	<b>DFIQKIANN</b>	<b>YACKHPEVQS</b>	<b>ILKISQPQEP</b>	<b>ELMNNANPSP</b>	300
301	<b>PSPSQINLG</b>	<b>PSSNPHAKPS</b>	<b>DFHFLKVIGK</b>	<b>GSFGKLLAR</b>	<b>HKAEEVFYAV</b>	<b>KVLQKKALK</b>	360
361	<b>KKEEKHIMSE</b>	<b>RNVLLKNVKH</b>	<b>PFLVGLHFSF</b>	<b>QTADKLYFVL</b>	<b>DYINGGELFY</b>	<b>HLQRERCFLE</b>	420
421	<b>PRARFYAAEI</b>	<b>ASALGYLHSL</b>	<b>NIVYRDLKPE</b>	<b>NILLDSQGHI</b>	<b>VLTDFGLCKE</b>	<b>NIEHNSTTST</b>	480
481	<b>FCGTPEYLPE</b>	<b>VLHKQPYDRT</b>	<b>VDWWCLGAVL</b>	<b>YEMLYGLPPF</b>	<b>YSRNTAEMD</b>	<b>NILNKPLQLK</b>	540
541	<b>PNITNSARHL</b>	<b>LEGLLQKDR</b>	<b>KRLGAKDDFM</b>	<b>EIKSHVFFSL</b>	<b>INWDDLINLK</b>	<b>ITPPFNPVNS</b>	600
601	<b>GPNDLRHFDP</b>	<b>EFTEEPVNS</b>	<b>IGKSPDSVLV</b>	<b>TASVKEAEEA</b>	<b>FLGFSYAPPT</b>	<b>DSFL</b>	660

1-218: GST **blue**: SGK1

SGK1 wt <sup>1</sup> Amino Acid Sequence							
1	<b>MTVKTEAAKG</b>	<b>TLTYSRMRGM</b>	<b>VAILIAFMKQ</b>	<b>RRMGLNDFIQ</b>	<b>KIANNYSYACK</b>	<b>HPEVQSILKI</b>	60
61	<b>SQPQEPPELMN</b>	<b>ANSPPPSPS</b>	<b>QQINLGSSN</b>	<b>PHAKPSDFHF</b>	<b>LKVIGKGSFG</b>	<b>KVLLARHKA</b>	120
121	<b>EVFYAVKVLQ</b>	<b>KKAILKKKEE</b>	<b>KHIMSERNL</b>	<b>LKNVKHPFLV</b>	<b>GLHFSFQTAD</b>	<b>KLYFVLDYIN</b>	180
181	<b>GGELFYHLQR</b>	<b>ERCFLEPRAR</b>	<b>FYAAEIASAL</b>	<b>GYLHSLNIVY</b>	<b>RDLKPENILL</b>	<b>DSQGHIVLTD</b>	240
241	<b>FGLCKENIEH</b>	<b>NSTTSTFCGT</b>	<b>PEYLAPEVLH</b>	<b>KQPYDRTVDW</b>	<b>WCLGAVLYEM</b>	<b>LYGLPPFYSR</b>	300
301	<b>NTAEMYDNIL</b>	<b>NKPLQLKPN</b>	<b>TNSARHLEG</b>	<b>LLQKDRTRKRL</b>	<b>GAKDDFMEIK</b>	<b>SHVFFSLINW</b>	360
361	<b>DDLINKKITP</b>	<b>PFNPNVSGPN</b>	<b>DLRHFDPPEFT</b>	<b>EEPVPNSIGK</b>	<b>SPDSVLVTAS</b>	<b>VKEAAEAFLG</b>	420
421	<b>FSYAPPTDSF</b>	<b>L</b>					480

**blue**: SGK1 sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](https://www.ncbi.nlm.nih.gov/protein/NP_005618.2) accession number NP\_005618.2