

## Certificate of Analysis

### SGK1 (S422D), unactive

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-332, 14-332-K, 14-332M

Parent Lot # 1792535

The data presented in this document apply to the parent lot shown above and to all pack sizes derived from subsequent vialling runs of this parent lot. An alphabetical suffix after the parent lot number is used to denote each vialling run.

**Product Description:** N-terminal 6His-tagged, recombinant, human SGK1, amino acids 60–end, containing the mutation S422D. Expressed by baculovirus in Sf21 insect cells. Purified using Ni<sup>2+</sup>/NTA agarose. Purity 97.3% by SDS-PAGE and Coomassie blue staining. MW = 45.7kDa.

**Specific Activity (Parent lot# 1792535):** 1451U/mg after activation with PDK1 as described below. One Unit of SGK1 (S422D) activity is defined as 1nmol phosphate incorporated into Crosstide in one minute.

**Formulation:** 2.935mg/ml of enzyme in 50mM Tris/HCl pH7.5, 0.1mM EGTA, 150mM NaCl, 0.03% Brij-35, 0.2mM PMSF, 1mM benzamidine, 0.1% 2-mercaptoethanol, 270mM sucrose. Frozen solution.

**Storage and Stability:** On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 6 months from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

**Handling Recommendations:** Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled microcentrifuge tubes and immediately snap-freeze the vials in liquid nitrogen prior to re-storage at -70°C

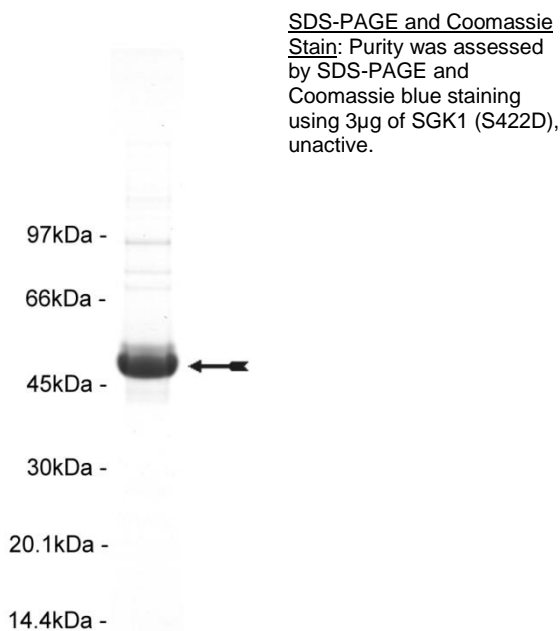
**FOR IN VITRO RESEARCH USE ONLY  
NOT FOR USE IN HUMANS OR ANIMALS**

### Quality Control Testing

**Activation Assay:** Unactive SGK1 (S422D) was activated using PDK1 (cat# 14-452) and the increased activity against crosstide determined. The activation and subsequent assay is described on page two. Results of this assay are shown below.

**MS Tryptic Fingerprint:** Confirmed identity as SGK1 (S422D) with the translated sequence listed on page three.

Active PDK1	Unactive SGK1 (S422D)	ng linear range	Mean cpm	Comments
10ng	500ng	5–20	7864 – 25505	Activated SGK1
2.5µg	7.4µg	50–200	340 – 716	Background



## Certificate of Analysis

### Kinase Cascade Assay Protocol

#### Stock Solutions:

1. **10 x SGK1 activation buffer:** 500mM Tris/HCl pH7.5, 1mM EGTA, 1% 2-mercaptoethanol.
2. **5 x SGK1 assay buffer:** 40mM MOPS/NaOH pH7.0, 1mM EDTA.
3. **Enzyme Dilution Buffer:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 5% glycerol, 0.01% Brij-35, 0.1% 2-mercaptoethanol, 1mg/ml BSA.
4. **Magnesium/ATP Cocktail (5 x stock):** 1mM cold ATP and 100mM magnesium acetate.
5. **[ $\gamma$ -<sup>33</sup>P]ATP:** 2.5 x magnesium acetate/[ $\gamma$ -<sup>33</sup>P]ATP cocktail: 25mM MgAc and 0.25mM ATP to which is added [ $\gamma$ -<sup>33</sup>P]ATP (specific activity approximately 500 - 800cpm/pmol as required.)
6. **PDK1, active (Catalogue# 14-452):** Use at a final assay concentration of 0.025 $\mu$ M. Prepare a 0.0147mg/ml stock and add 2.5 $\mu$ l of stock per assay point.
7. **SGK1 (S422D), unactive:** Use at a final concentration of 4 $\mu$ M (0.1828mg/ml). Prepare a 0.914mg/ml stock and add 5 $\mu$ l of stock per assay point.
8. **Modified Crosstide (GRPRTSSFAEGKK):** Use a final assay concentration of 30 $\mu$ M. Make a 300 $\mu$ M stock. Add 2.5 $\mu$ l of stock per assay point.

#### Assay Procedure:

##### **Stage One:** *Activation of SGK1 (S422D) by PDK1.*

1. Add 5 $\mu$ l of SGK1 activation buffer to a microcentrifuge tube.
2. Add 5 $\mu$ l of **PDK1, active.**
3. Add **5 $\mu$ l of SGK1 (S422D), unactive.**
4. Add 30 $\mu$ l of dH<sub>2</sub>O.
5. Add 5 $\mu$ l of magnesium/ATP cocktail.
6. Incubate for 30 minutes at 30°C.
7. Stop reaction by diluting 10–50 fold and placing on ice.

##### **Stage Two:** *Phosphorylation of modified crosstide by SGK1 (S422D) (96 well plate format)*

1. Add 5 $\mu$ l of 5 x reaction buffer per assay to wells
2. Add 2.5 $\mu$ l of **modified Crosstide (GRPRTSSFAEGKK).**
3. Add **2.5 $\mu$ l (5–20ng) SGK1 (S422D), active from Stage One.**
4. Add 5 $\mu$ l of dH<sub>2</sub>O.
5. Add 10 $\mu$ l of diluted [ $\gamma$ -<sup>33</sup>P] ATP mixture.
6. Incubate for 10 minutes at 30°C.
7. Stop the reaction by adding 5 $\mu$ l of 3% phosphoric acid.
8. Transfer a 10 $\mu$ l aliquot onto the appropriate area of a **P30 Filtermat.**
9. Wash the filtermat twice for 5 minutes with 75mM phosphoric acid.
10. Wash the filtermat once for 2 minutes with methanol.
11. Transfer the filtermat to a sealable plastic bag and add 4ml of scintillation cocktail.
12. Read in a scintillation counter. Compare cpm of enzyme samples with cpm of control samples that contain all assay components plus 1 $\mu$ l of 30% phosphoric acid.

## Certificate of Analysis

### SGK1 (S422D) Sequence Information

<b><u>Protein</u></b>	Human SGK1 (S422D)
<b><u>Tags</u></b>	N-terminal 6His
<b><u>Native sequence</u></b>	I28 of the recombinant protein is equivalent to I60 of human SGK1
<b><u>Accession number</u></b>	GenBank AF153609

#### Recombinant SGK1 (S422D) amino acid sequence:

```

1 MSYYHHHHHH DYDIPTTENL YFQGAMGISQ PQEPELMNAN PSPPPSPSQQ INLG PSSNPH
61 AKPSDFHFLK VIGKGSFGKV LLARHKAEEV FYAVKVLQKK AILKKKEEKH IMSERNVLLK
121 NVKHPFLVGL HFSFQTADKL YFVLDYINGG ELFYHLQRRER CFLEPRARFY AAEIASALGY
181 LHSLNIVYRD LKPENILLDS QGHIVLTDG LCKENIEHNS TTSTFCGTP EYL APEVLHKQ
241 PYDRTVDWWC LGAVLYEMLY GLPPFYSRNT AEMYDNILNK PLQLKPNITN SARHLLLEGLL
301 QKDRTKRLGA KDDFMEIKSH VFFSLINWDD LINKKITPPF NPNVSGPNDL RHFDP EFTEE
361 PVPNSIGKSP DSVLVTASVK EAAEAF LGFD YAPPTDSFL
  
```

#### Recombinant SGK1 (S422D) nucleotide sequence:

```

1 atgtcgtact accatcacca tcaccatcac gattacgata tcccaacgac cgaaaacctg
61 tattttcagg ggcctatggg gatctcccaa cctcaggagc ctgagcttat gaatgccaac
121 ccttctcctc caccaagtcc ttctcagcaa atcaaccttg gcccgctgct caatcctcat
181 gctaaacctc ctgactttca cttcttgaaa gtgatcggaa agggcagttt tggaaagggt
241 cttctagcaa gacacaaggc agaagaagtg ttctatgcag tcaaagtttt acagaagaaa
301 gcaatcctga aaaagaaaga ggagaagcat attatgtcgg agcggaatgt tctgttgaag
361 aatgtgaagc accctttcct ggtgggcctt cacttctctt tccagactgc tgacaaattg
421 tactttgtcc tagactacat taatggtgga gagttgttct accatctcca gagggaacgc
481 tgcttcctgg aaccacgggc tcgtttctat gctgctgaaa tagccagtgc cttgggctac
541 ctgcattcac tgaacatcgt ttatagagac ttaaaccagc agaatatttt gctagattca
601 cagggacaca ttgtccttac tgatttcgga ctctgcaagg agaacattga acacaacagc
661 acaacatcca cttctgtggt cacgccggag tatctcgcac ctgaggtgct tcataagcag
721 cttatgaca ggactgtgga ctgggtggtgc ctgggagctg tcttgatga gatgctgat
781 ggctgcccgc ctttttatag ccgaaacaca gctgaaatgt acgacaacat tctgaacaag
841 cctctccagc tgaacacaaa tattacaaat tccgcaagac acctcctgga gggcctcctg
901 cagaaggaca ggacaaagcg gctcggggcc aaggatgact tcatggagat taagagtcac
961 gtcttcttct ccttaattaa ctgggatgat ctcattaata agaagattac tccccctttt
1021 aacccaaatg tgagtgggac caacgacctc cggcactttg accccgagtt taccgaagag
1081 cctgtcccca actccattgg caagtcccct gacagcgtcc tcgtcacagc cagcgtcaag
1141 gaagctgccg aggtttcctt aggtttgac tatgcgcctc ccacggactc tttcctctga
  
```

Reviewed and approved by site quality representative.

Unless otherwise stated in our catalogue or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

© 2014 Eurofins Pharma Discovery Services UK Limited is an independent member of Eurofins Discovery Services.