

## ProQinase™ TSSK3

testis specific serine kinase 3

### Recombinant Human Active Protein Kinase

HGNC Symbol: TSSK3

Synonyms: STK22C, SPOGA3

Product No.: 2060-0000-1

Lot: 004

**Description:** Human TSSK3, full length, amino acids M<sub>1</sub>-T<sub>268</sub> (as in [NCBI/Protein](#) entry NP\_443073.1), untagged, expressed in Sf9 insect cells

**Product identity:** TSSK3 Lot 004, has been verified by mass spectrometry LC-ESI-MS/MS

**Theoretical MW**<sub>Fusion Protein</sub>: 30256 Da

**Expression host:** Sf9 insect cells

**Purification:** GST-Affinity Chromatography

**Activation:** This kinase was not activated by special procedures

**Storage buffer:** 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 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!

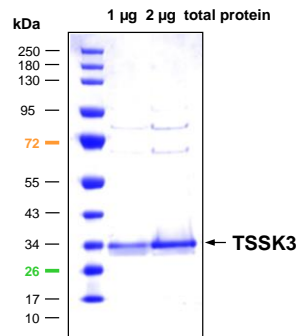
**Protein concentration:** 0.091 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Biochemical Parameters:**

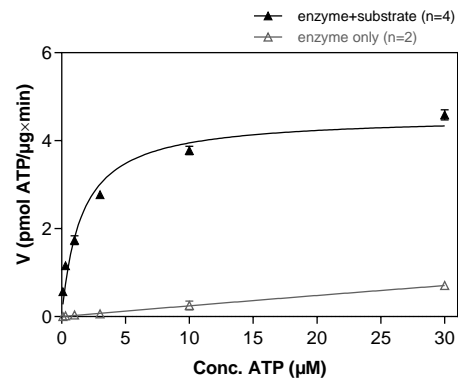
Specific kinase activity (P<sub>i</sub> transfer): 4.6 pmol/µg\*min

ATP-K<sub>M</sub>: 1.5 µM

**TSSK3 Lot 004:  
Coomassie stain**



**TSSK3 Lot 004:  
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: 4xChocktide 10 µg/ml
  - Kinase: 6 µg/ml

Assay technology:  
Radiometric filter binding assay  
MSFC membrane (96 well plate, Millipore)

Recombinant Proteins

Sequence information

GST-TSSK3 Recombinant Fusion Protein Amino Acid Sequence								
1	GP	MEDFLLSN	GYQLGKTIGE	GTYSKVKEAF	SKKHQRKVAI	KVIDKMGGPE	EFIQRFLPRE	60
61	LQ	IVRTL DHK	NIIQVYEMLE	SADGKICLVM	ELAEGGDVFD	CVLNGGPLPE	SRAKALFROM	120
121	VE	AIRYCHGC	GVAHRDLKCE	NALLQGFLNK	LTDFGFAKVL	PKSHRELSQT	FCGSTAYAAP	180
181	EV	LQGI PHDS	KKGDVWSMGV	VLYVMLCASL	PFDDTDIPKM	LWQQQKGVSF	PTHLSISADC	240
241	QD	LLKRLLEP	DMILRPSIEE	VSWHPWLAST				300

1-2: legacy of 3C protease cleavage **Red:** HIS6-tag **blue:** TSSK3

TSSK3 wt <sup>1</sup> Amino Acid Sequence								
1	MED	FLLSNGY	QLGKTIGEGT	YSKVKEAFSK	KHQRKVAIKV	IDKMGGPPEF	IQRFLPRELQ	60
61	IVR	TL D HKNI	IQVYEMLESA	DGKICLVMEL	AEGGDVFD CV	LNGGPLPESR	AKALFROMVE	120
121	AIR	YCHGCGV	AHRDLKCENA	LLOGFNLKL T	DFGFAKVL PK	SHRELSQTFC	GSTAYA APEV	180
181	LQ	GI PHDSK	KGDVWSMGVVL	YVMLCASL PF	DDTDIPKMLW	QQQKGVSFPT	HLSISADCQD	240
241	LL	KRLLEPDM	ILRPSIEEVS	WHPWLAST				300

**blue:** TSSK3 sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](https://www.ncbi.nlm.nih.gov/protein/NP_443073.1) accession number NP\_443073.1