

ProQinase™ TSF1

serine/threonine kinase 16

Recombinant Human Active Protein Kinase

HGNC Symbol: STK16

Synonyms: FLJ39635, hPSK, KRCT, MGC16211, MPSK, MPSK1, PKL12

Product No.: 0187-0000-1

Lot: 002

Description: Human TSF1, full length, amino acids M₁-I₃₀₅ (as in [NCBI/Protein](#) entry NP_001008910.1), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: TSF1 Lot 002, was confirmed as TSF1 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 64,393 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: This kinase was not activated by special procedures

Storage buffer: 50 mM TRIS-HCl pH 8.0, 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: 1.02 µg/µl

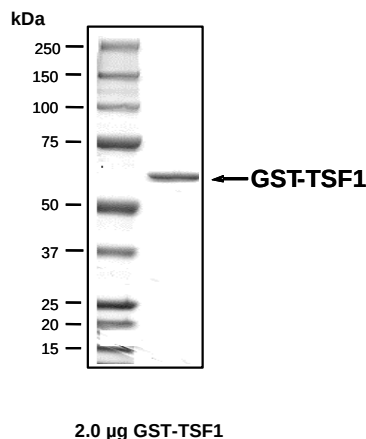
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

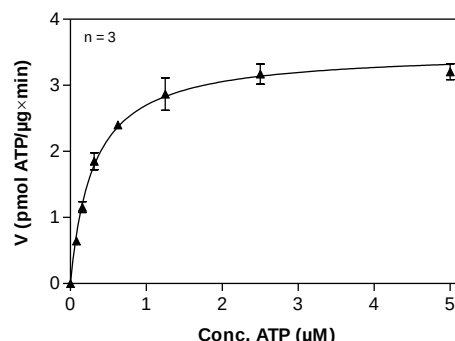
Specific kinase activity (P_i transfer): 3.5 pmol/µg × min

ATP-K_M: 0.3 µM

TSF1 Lot 002: Coomassie stain



TSF1 Lot 002: 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: Casein 200 µg/ml
Kinase: 2 µg/ml
- Filter binding assay
MSFC membrane (Millipore)

Additional assay technology:

TSF1 Lot 002 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

ProQinase™ TSF1

Product No.: 0187-0000-1

GST-TSF1 Recombinant Fusion Protein Amino Acid Sequence

1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLIERDEGDK	WRNKKFELGL	EFNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHH	RRRASVAAGI	240
241	LVPRGSPGLD	GIYARGIQMG	HALCVCSSGT	VIIDNKRYLF	IQKLGEFFS	YVDLVEGLHD	300
301	GHFYALKRIL	CHEQQDREEA	QREADMHRLE	NHPNLRILVA	YCLRERGAH	EAWLLLPFFK	360
361	RGTWNEIER	LKDKGNFLE	DQILWLLGI	CRGLEAIHAK	GYAHRDLKPT	NILLGDEGQP	420
421	VLMDLGSNNQ	ACIHVEGSRQ	ALTLDWAAQ	RCTISYRAPE	LFSVQSHCVI	DETDVWSLG	480
481	CVLYAMMFGE	GPYDMVFQKG	DSVALAVQNG	LSIPQSPRHS	SALRQLNSM	MTVDPHQRPH	540
541	IPLLLSQLEA	LQPPAPGQHT	TQI				600

1-218: GST **Red**: HIS6-tag **Pink**: Thrombin cleavage site **blue**: TSF1

TSF1 wt¹ Amino Acid Sequence

1	MGHALCVCSSR	GTVIIDNKRY	LFIQKLGEFG	FSYVDLVEGL	HDGHFYALKR	ILCHEQQDRE	60
61	EAQREADMHR	LFNHPNLRIL	VAYCLRERGA	KHEAWLLLPF	FKRGTWNEI	ERLKDKNFL	120
121	TEDQILWLLI	GICRGLEATH	AKGYAHRDLK	PTNILLGDEG	QPVLMDLGSM	NQACIHVEGS	180
181	RQALTLDWA	AQRCTISYRA	PELFSVQSHC	VIDERTDVWS	LGCVLYAMMF	GEGPYDMVFQ	240
241	KGDSVALAVQ	NQLSIPQSPR	HSSALRQLLN	SMMTVDPHQR	PHIPLLLSQL	EALQPPAPGQ	300
301	HTTQI						360

blue: TSF1 sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_001008910.1