

## ProQinase™ MST1

serine/threonine kinase 4 (STK4)

Recombinant Human Active Protein Kinase

HGNC Symbol: STK4

Synonyms: KRS2, YSK3

Product No.: 0667-0000-1

Lot: 002

**Description:** Human MST1, amino acids T<sub>3</sub>-F<sub>487</sub> (as in [NCBI/Protein](#) entry NP\_006273.1), N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

**Product identity:** MST1, Lot 002, was confirmed as MST1 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW**<sub>Fusion Protein</sub>: 86,340 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, 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:** 0.306 µg/µl

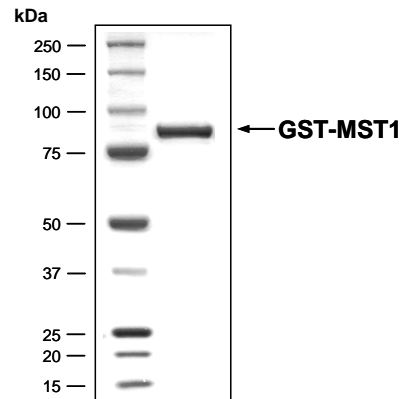
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Biochemical Parameters:**

Specific kinase activity (P<sub>i</sub> transfer): 322 pmol/µg × min

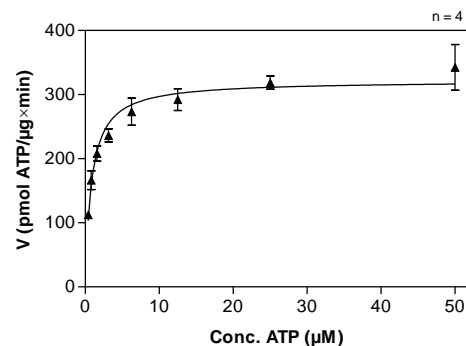
ATP-K<sub>M</sub>: 0.8 µM

### MST1 Lot 002: Coomassie stain



2.0 µg GST-MST1

### MST1 Lot 002: Determination of V<sub>max</sub> and K<sub>M</sub> value for ATP



### Determination of K<sub>M</sub> value & Specific activity:

- 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: Myelin Basic Protein 100 ng/ml
- MST1: 0.5 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

### Additional assay technology:

MST1 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™ MST1

Product No.: 0667-0000-1

GST-MST1 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	PMGHHHHHG	RRRASVAAGI	240
241	LVPRGSPGLD	GIYARDSLEV	LFQGPLAMLT	VQLRNPPRRQ	LKKLDEDSL	KQPEEVFDVL	300
301	EKLGEFSYGS	VYKAIHKETG	QIVAIKQVPV	ESDLQEIIKE	ISIMQQCDSP	HVVKYGSYF	360
361	KNTDLWIVME	YCGAGSVSDI	IRLRNKTLTE	DEIATILQST	LKGLEYLHFM	RKIHRDIKAG	420
421	NILLNTEGHA	KLADFGVAGQ	LTDMAKRNT	VIGTPFWMAP	EVIQEIGYNC	VADIWSLGIT	480
481	AIEMAEGKPP	YADIHPMRAI	FMIPTNPPPT	FRKPELWSDN	FTDFVKQCLV	KSPEQRATAT	540
541	QLLQHPFVRS	AKGVSILRDL	INEAMDVKLK	RQESQREVD	QDDEENSEED	EMDSGTMVRA	600
600	VGDEMGTVRV	ASTMTDGANT	MIEHDDTLPS	QLGTMVINA	DEEEEGTMKR	RDETMQPAK	660
661	SFLEYFEQHE	KENQINSFGK	SVPGPLKNS	DWKIPQGDY	EFLKSWTVED	LQKRLALDP	720
721	MMEQEIEEIR	QKYQSKRQPI	LDAIEAKRR	QQNF			780

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site Green: 3C cleavage site blue: MST1 boxed: variation from RefSeq

MST1 wt <sup>1</sup> Amino Acid Sequence							
1	METVQLRNPP	RRQLKKLDED	SLTKQPEEVF	DVLEKLGEKS	YGSVYKAIHK	ETGQIVAIKQ	60
61	VPVESDLQEI	IKEISIMQQC	DSPHVVKYYG	SYFKNTDLWI	VMEYCGAGSV	SDIIRLRNKT	120
121	LTEDEIATIL	QSTLKGLEYL	HFMRKIHRDI	KAGNILLNTE	GHAKLADFGV	AGQLTDTMAK	180
181	RNTVIGTPFW	MAPEVIQEIG	YNCVADIWSL	GITAIEMAEG	KPPYADIHPM	RAIFMIPTNP	240
241	PPTFRKPELW	SDNFTDFVKQ	CLVKSPEQRA	TATQLLQHPF	VRSAGVSIL	RDLINEAMDV	300
301	KLKRQESQQR	EMDQDDEENS	EEDEMSGTM	VRAVGDEMGT	VRVASTMTDG	ANTMIEHDDT	360
361	LPSQLGTMVI	NAEDEEEEGT	MKRRDETMQP	AKPSFLEYFE	QKEKENQINS	FGKSVGPLK	420
421	NSSDWKIPQD	GDYEFKLSWT	VEDLQKRLLA	LDPMMEQEIE	EIRQKYQSKR	QPILDAIEAK	480
481	KRRQNF						540

blue: MST1 sequence expressed in recombinant protein Red: variant in recombinant protein

<sup>1</sup>NCBI/Protein accession number NP\_006273.1

M312V: SNP variation see NCBI/dbSNP ID: rs17420378