

## MST2

serine/threonine kinase 3

Recombinant Human Active Protein Kinase

HGNC Symbol: STK3

Synonyms: KRS1

Product No.: 0699-0000-1

Lot: 003

**Description:** Human MST2, full length, amino acids M<sub>1</sub>-F<sub>491</sub> (as in [NCBI/Protein](#) entry NP\_006272.2), untagged, expressed in Sf9 insect cells

**Product identity:** MST2 Lot 003, was confirmed as MST2 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW**<sub>Fusion Protein</sub>: 56,884 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:** 0.271 µg/µl

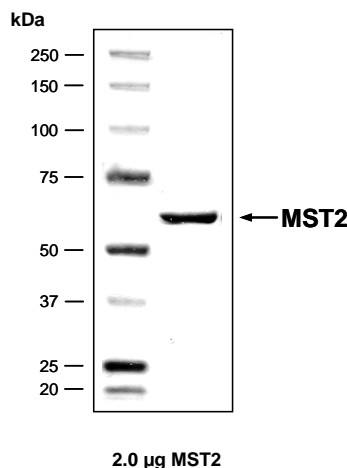
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Biochemical Parameters:**

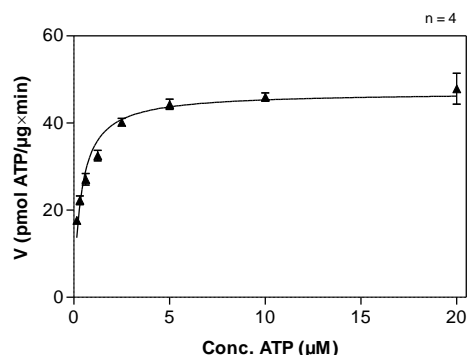
Specific kinase activity (P<sub>i</sub> transfer): 47 pmol/µg × min

ATP-K<sub>M</sub>: 0.38 µM

**MST2 Lot 003:  
Coomassie stain**



**MST2 Lot 003:  
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: Casein 100 µg/ml
  - Kinase: 1 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

**Additional assay technology:**

MST2 Lot 003 was also successfully tested by ProQinase 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

This product was manufactured at ProQinase in Freiburg, Germany, and is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.

## MST2

Product No.: 0699-0000-1

MST2 Recombinant Fusion Protein Amino Acid Sequence								
1	GPLAML	MEQP	PAPKSKLKKL	SEDSLTKQPE	EVFDVLEKLG	EGSYGSVFKA	IHKESGQVVA	60
61	IKQVPVESDL	QEIIKEISIM	QQCDSPYVVK	YYGSYFKNTD	LWIVMEYCGA	GSVSDIIRLR		120
121	NKTLIEDEIA	TILKSTLKGL	EYLHFMRKIH	RDIKAGNILL	NTEGHAKLAD	FGVAGQLTDT		180
181	MAKRNTVIGT	PFWMAPEVIQ	EIGYNCVADI	WSLGITSIEM	AEGKPPYADI	HPMRAIFMIP		240
241	TNPPPTFRKP	ELWSDDFTFD	VKKCLVKNPE	QRATATQLLQ	HPFIKNAKPV	SILRDLITEA		300
301	MEIKAKRHEE	QQRELEEEEE	NSDEDELDSH	TMVKTSVESV	GTMRATSTMS	EGAQTMIEHN		360
361	STMLESDLGT	MVINSEDEEE	EDGTMKRNAT	SPQVQRPSFM	DYFDKQDFKN	KSHENCNQNM		420
421	HEPFPMSKNV	FPDNWKVPQD	GDFDFLKNLS	LEELQMLKA	LDPMMEREIE	ELRQRYTAKR		480
481	QPILDAMDAK	KRRQQNF						540

1-6: legacy of tag cleavage blue: MST2

MST2 wt <sup>1</sup> Amino Acid Sequence							
1	MEQPPAPKSK	LKKLSEDSL	KQPEEVFDVL	EKLGEESYGS	VFKAIHKESG	QVVAIKQVPV	60
61	ESDLQEIIKE	ISIMQQCDSP	YVVKYGSYF	KNTDLWIVME	YCGAGSVSDI	IRLRNKTLIE	120
121	DEIATILKST	LKGLEYLHFM	RKIHRDIKAG	NILLNTEGHA	KLADFGVAGQ	LTDITMAKRNT	180
181	VIGTPFWMAP	EVIQEIGYNC	VADIWLSGIT	SIEMAEGKPP	YADIHPMRAI	FMIPTNPPPT	240
241	FRKPELWSD	FTDFVKKCLV	KNPEQRATAT	QLLQHPFIKN	AKPVSILRDL	ITEAMEIKAK	300
301	RHEEQQRELE	EEEEENDEDE	LDSTMVKTS	VESVGTMRAT	STMSEGAQTM	IEHNSTMLES	360
361	DLGTMVINSE	DEEEEDGTMK	RNATSPQVQR	PSFMDYFDKQ	DFKNKSHENC	NQNMHEPFPM	420
421	SKNVFPDNWK	VPQDGFDFL	KNLSLEELQM	RLKALDPMME	REIEELRQRY	TAKRQPILDA	480
481	MDAKRRQQN	F					540

blue: MST2 sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](#) accession number NP\_006272.2

This product was manufactured at ProQinase in Freiburg, Germany, and is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.