

ProQinase™ 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 M₁-F₄₉₁ (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_{Fusion Protein}: 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

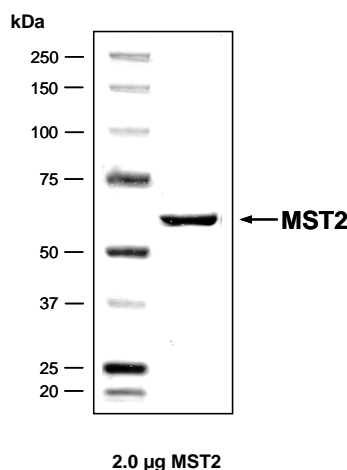
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

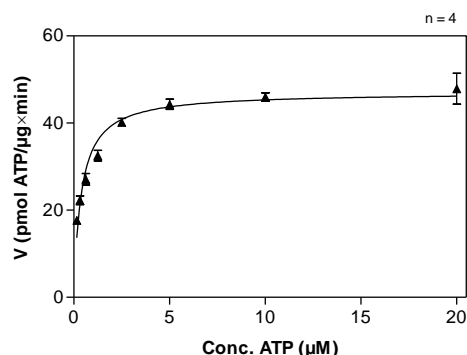
Specific kinase activity (P_i transfer): 47 pmol/µg × min

ATP-K_M: 0.38 µM

**MST2 Lot 003:
Coomassie stain**



**MST2 Lot 003:
Determination of V_{max} and K_M value for ATP**



Determination of K_M value & Specific activity:

- 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 100 µg/ml
 - Kinase: 1 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

Additional assay technology:

MST2 Lot 003 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™ MST2

Product No.: 0699-0000-1

MST2 Recombinant Fusion Protein Amino Acid Sequence								
1	GPLA	MEQP	PAPKSKLKKL	SEDSLTKQPE	EVFDVLEKLG	EGSYGSVFKA	IHKESGQVVA	60
61	IKQVP	VESDL	QEIIKEISIM	QQCDSPYVVK	YYGSYFKNTD	LWIVMEYCGA	GSVSDIIRLR	120
121	NKTLIE	DEIA	TILKSTLKGL	EYLHFMRKIH	RDIKAGNILL	NTEGHAKLAD	FGVAGQLTDT	180
181	MAKRNT	VIGT	PFWMAPEVIQ	EIGYNCVADI	WSLGITSIEM	AEGKPPYADI	HPMRAIFMIP	240
241	TNPPPT	FRKP	ELWSDDFTFD	VKKCLVKNPE	QRATATQLLQ	HPFIKNAKPV	SILRDLITEA	300
301	MEIKAKR	HEE	QQRELEEEEE	NSDEDELDSH	TMVKTSVESV	GTMRATSTMS	EGAQTMIEHN	360
361	STMLES	DLGT	MVINSEDEEE	EDGTMKRNAT	SPQVQRPSFM	DYFDKQDFKN	KSHENCNQNM	420
421	HEPFPMS	KNV	FPDNWKVPQD	GDFDFLKNLS	LEELQMLKA	LDPMMEREIE	ELRQRYTAKR	480
481	QPILDAM	DAK	KRRQQNF					540

1-6: legacy of tag cleavage blue: MST2

MST2 wt ¹ Amino Acid Sequence								
1	MEQP	PAPKSK	LKKLSE	DSLTKQPEE	VFDVLEKLG	EGSYGSVFKA	IHKESGQVVA	60
61	ESDL	QEIIKE	ISIMQQ	CDSPYVVK	YYGSYFKNTD	LWIVMEYCGA	GSVSDIIRLR	120
121	DEIA	TILKST	LKGLYLFHM	RKIHRDIKAG	NILLNTEGHA	KLADFGVAGQ	LTDITMAKRNT	180
181	VIGT	PFWMAP	EVIQEIGYNC	VADIWLSGIT	SIEMAEGKPP	YADIHPMRAI	FMIPTNPPPT	240
241	FRKPEL	WSD	FTDFVKKCLV	KNPEQRATAT	QLLQHPFIKN	AKPVSILRDL	ITEAMEIKAK	300
301	RHEEQ	QRELE	EEEE	NSDEDE	LDSTMVKTS	VESVGTMRAT	STMSEGAQTM	360
361	DLGTM	VINSE	DEEEDG	TMKRNAT	SPQVQR	PSFMDYFDKQ	DFKNKSHENC	420
421	SKNV	FPDN	WKVPQD	GDFDFL	KNLSLEELQ	MLKALDPMME	REIEELRQRY	480
481	MDAK	KRRQQ	NF					540

blue: MST2 sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_006272.2