

ProQinase™ MLK3

mitogen-activated protein kinase kinase kinase 11

Recombinant Human Active Protein Kinase

HGNC Symbol: MAP3K11

Synonyms: PTK1, SPRK, MEKK11

Product No.: 1999-0000-1

Lot: 001

Description: Human MLK3, N-terminal fragment, amino acids M₁-G₄₈₈ (as in [NCBI/Protein](#) entry NP_002410.1), N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: MLK3 Lot 001, has been verified by mass spectrometry LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 82377 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, 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!

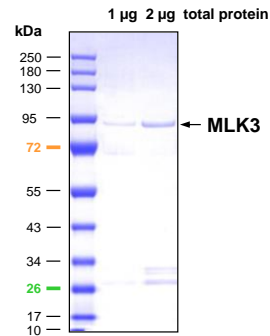
Protein concentration: 0.52 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

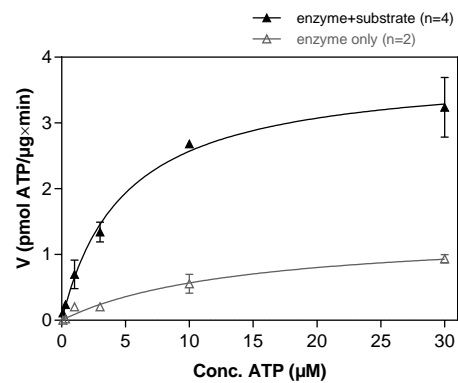
Specific kinase activity (P_i transfer): 3.8 pmol/µg*min

ATP-K_M: 4.9 µM

MLK3 Lot 001: Coomassie stain



MLK3 Lot 001: 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: Myelin basic protein 20 µg/ml
- Kinase: 1 µg/ml

Assay technology:

- Radiometric filter binding assay
- MSFC membrane (96 well plate, Millipore)

Recombinant Proteins

Sequence information

| GST-MLK3 Recombinant Fusion Protein Amino Acid Sequence | | | | | | | |
|---|-------------|------------|------------|------------|------------|------------|-----|
| 1 | MSPILGYWKI | KGLVQPTRL | LEYLEEKYEE | HLYERDEGDK | WRNKKFELGL | EFPNLPYYID | 60 |
| 61 | GDVKLTQSMA | IIRYIADKHN | MLGGCPKERA | EISMLEGAVL | DIRYGVSRIA | YSKDFETLKV | 120 |
| 121 | DFLSKLPPEML | KMFKDRLCHK | TYLNGDHVTH | PDFMLYDALD | VVLYMDPMCL | DAFPKLVCFK | 180 |
| 181 | KRIEAIPIQID | KYLKSSKYIA | WPLQGWQATF | GGGDHPPKSD | PMGHHHHHGG | RDSLEVLFGG | 240 |
| 241 | MEPLKSLFL | KSPLGSWNGS | SGGGGGGGGG | GRPEGSPKAA | GYANPVWTAL | FDYEPSGQDE | 300 |
| 301 | LALRKGDRVE | VLSRDAAISG | DEGWAGQVG | QVGIFFPSNY | VSRGGGPPPC | EVASFQELRL | 360 |
| 361 | EEVIGIGGFG | KVYRGSWRGE | LVAVKAARQD | PDEDISVTAE | SVRQEARLFA | MLAHPNIIAL | 420 |
| 421 | KAVCLEEPNL | CLVMEYAAGG | PLSRALAGR | VPPHVLVNW | VQIARGMHYL | HCEALVPVIH | 480 |
| 481 | RDLKSNNILL | LQPIESDDME | HKTLKITDFG | LAREWHKTQ | MSAAGTYAWM | APEVIKASTF | 540 |
| 541 | SKGSDVWSFG | VLLWELLTGE | VPYRGIDCLA | VAYGVAVNKL | TLPIPSTCPE | PFAQLMADCW | 600 |
| 601 | AQDPHRRPDF | ASILQOLEAL | EAQVLEMPR | DSFHSMQEGW | KREIQGLFDE | LRAKEKELLS | 660 |
| 661 | REEELTRAAR | EQRSQAEQLR | RREHLLAQWE | LEVFERELTL | LLQQVDRERP | HVRRRRGTFK | 720 |
| 721 | RSKLARDG | | | | | | 780 |

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: MLK3 fragment

| MLK3 wt ¹ Amino Acid Sequence | | | | | | | |
|--|------------|-------------|-------------|--------------|------------|------------|-----|
| 1 | MEPLKSLFLK | SPLGSWNGSG | SGGGGGGGGG | RPEGSPKAAG | YANPVWTALF | DYEPSGQDEL | 60 |
| 61 | ALRKGDRVEV | LSRDAAISGD | EGWWAGQVGG | QVGIFFPSNYV | SRGGGPPPC | VASFQELRLE | 120 |
| 121 | EVIGIGGFGK | VYRGSWRGEL | VAVKAARQDP | DEDISVTAES | VRQEARLFAM | LAHPNIIALK | 180 |
| 181 | AVCLEEPNLC | LVMEYAAGGP | LSRALAGR | VPPHVLVNWAV | QIARGMHYH | CEALVPVIHR | 240 |
| 241 | DLKSNNILL | LQPIESDDMEH | KTLKITDFGL | AREWHKTQ | SAAGTYAWMA | PEVIKASTFS | 300 |
| 301 | KGSDVWSFGV | LLWELLTGEV | PYRGIDCLAV | AYGVAVNKL | LPIPSTCPEP | FAQLMADCWA | 360 |
| 361 | QDPHRRPDFA | SILQOLEALE | AQVLEMPRD | SFHSMQEGWK | REIQGLFDEL | RAKEKELLSR | 420 |
| 421 | EEELTRAARE | QRSQAEQLRR | REHLLAQWEL | EVFERELTLL | LQQVDRERPH | VRRRRGTFKR | 480 |
| 481 | SKLRARDGGE | RISMPDFKH | RITVQASPL | DRRRNVFEVG | PGDSPTFPRF | RAIQLEPAEP | 540 |
| 541 | GQAWGRQSPR | RLEDSSNGER | RACWAWGPSS | PKPGEAQNGR | RRSRMDEATW | YLDSDSSPL | 600 |
| 601 | GSPSTPPALN | GNPPRPSLEP | EEPKRVPVPAE | RGSSSGTPKL | IQRALLRGTA | LLASLGLGRD | 660 |
| 661 | LQPPGGPGRE | RGESPTTPPT | PTPAPCPTEP | PPSPLICFSL | KTPDSPPTPA | PLLLDLGIPV | 720 |
| 721 | GQRSASPRR | EEPRGGTVS | PPPGTSR | SAPGTPGTPRSP | LGLISRPRPS | PLRSRIDPWS | 780 |
| 781 | FVSAGPRPSP | LPSPQPAPRR | APWTLFPDSD | PFWDSPANP | FQGGPQDCRA | QTKDMGAQAP | 840 |
| 841 | WVPEAGP | | | | | | 900 |

blue: MLK3 sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_002410.1