

ProQinase™ MARK3

MAP/microtubule affinity-regulating kinase 3

Recombinant Human Active Protein Kinase

HGNC Symbol: MARK3

Synonyms: cTAK1, EMK2, KP78, PAR1A

Product No.: 0433-0000-1

Lot: 001

Description: MARK3, full length, amino acids M₁-L₇₁₃ (as in NCBI/Protein entry NP_001122392.2), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: MARK1 Lot 001, was confirmed as MARK1 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 109,443 Da

Expression: Baculovirus infected Sf9 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: 0.082 µg/µl

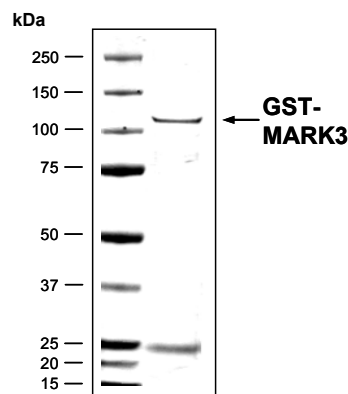
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

Specific kinase activity (P_i transfer): 18 pmol/µg × min

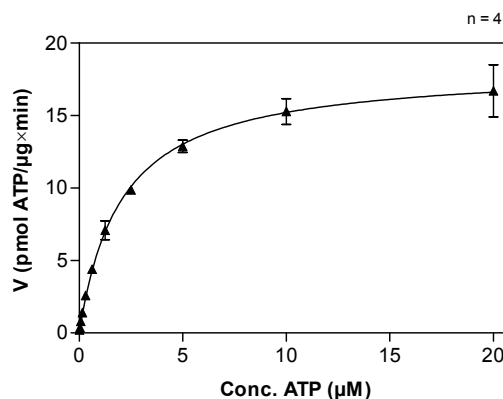
ATP-K_M: 2 µM

MARK3 Lot 001: Coomassie stain



2.0 µg GST-MARK3

MARK3 Lot 001: 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: Example#1 0.0 µg/ml
 - Kinase: 1.0 µg/ml
- Filter binding assay
 - MSFC/PH membrane (Millipore)

Additional assay technology: MARK3 Lot 001

was also successfully tested by Reaction Biology for the use with the ADP-Glo™ Kinase assay from ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details



ProQinase™ MARK3

Product No.: 0433-0000-1

| MARK3 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 | GIYARQMSTR | TPLPTVNERD | TENHTSHGDG | RQEVTSRTSR | SGARCRNSIA | 300 |
| 301 | SCADEQPHIG | NYRLKLTIGK | GNFAKVKLAR | HILTGREVAI | KIIDKTQLNP | TSLQKLFREV | 360 |
| 361 | RIMKILNHPN | IVKLFVET | EKTLYLIMEY | ASGGEVFDYL | VAHGRMKEKE | ARSKFRQIVS | 420 |
| 421 | AVQYCHQKRI | VHRDLKAENL | LLDADMNKI | ADFGFSNEFT | VGGKLDTFCG | SPPYAAPELF | 480 |
| 481 | QGKKYDGPEV | DVWSLGVILY | TLVSGSLPFD | GQNLKELRER | VLRGKYRIPF | YMSTDCENLL | 540 |
| 541 | KRFLVLNPIK | RGTLQIMKD | RWINAGHEED | ELKPFVEPEL | DISDQKRIDI | MVGMGYSQEE | 600 |
| 601 | IQESLSKMKY | DEITATYLLL | GRKSSEVRPS | SDLNNSGQS | PHHKVQRSVS | SSQKQRRYS | 660 |
| 661 | HAGPAIPSVV | AYPKRSQTST | ADSDLKEDGI | SSRKSSGSAV | GGKGIAPASP | MLGNASPNK | 720 |
| 721 | ADIPERKKSS | TVPSSNTASG | GMTRRNTYVC | SERTTADRHS | VIQNGKENST | IPDQRTPVAS | 780 |
| 781 | THSISSAATP | DRIRFPRGTA | SRSTFHGQPR | ERRTATYNGP | PASPPLSHEA | TPLSQTRSRG | 840 |
| 841 | STNLFSLKTS | KLTRSRNVSA | EQKDENEKAK | PRSLRFTWSM | KTTSSMDPGD | MMREIRKVL | 900 |
| 901 | ANNCDYEQRE | RELLFCVHGD | GHAENLVQWE | MEVCKLPRLS | LNGVRFKRIS | GTSIAFKNIA | 960 |
| 961 | SKIANELK | | | | | | 1020 |

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site blue:MARK3

| MARK3 wt ¹ Amino Acid Sequence | | | | | | | |
|---|-------------|------------|------------|------------|------------|------------|-----|
| 1 | MSTRTPPLPTV | NERDTENHTS | HGDGRQEVTS | RTSRSGARCR | NSIASCADEQ | PHIGNYRLK | 60 |
| 61 | TIGKGNFAKV | KLARHILTGR | EVAIKIIDKT | QLNPTSLQKL | FREVRIMKIL | NHPNIVKLFE | 120 |
| 121 | VIETEKTLYL | IMEYASGGEV | FDYLVAHGRM | KEKEARSKFR | QIVSAVQYCH | QKRIVHRDLK | 180 |
| 181 | AENLLLDADM | NIKIADFGFS | NEFTVGGKLD | TFCGSPPYAA | PELFQGGKYD | GPEVDVWSLG | 240 |
| 241 | VILYTLVSGS | LPFDGQNLKE | LRERVLRGKY | RIPFYMSTDC | ENLLKRFLVL | NPIKRGTLQ | 300 |
| 301 | IMKDRWINAG | HEEDELKPFV | EPELDIDSDQ | RIDIMVGMGY | SQEEIQESLS | KMKYDEITAT | 360 |
| 361 | YLLLGKRSSE | VRPSSDLNNS | TGQSPHKKVQ | RSVSSSQQR | RYSDHAGPAI | PSVVAYPKRS | 420 |
| 421 | QTSTADSDLK | EDGISSRKSS | GSAVGGKZIA | PASPMLGNAS | NPNKADIPER | KKSSTVPSSN | 480 |
| 481 | TASGGMTRRN | TYVCSERTTA | DRHSVIQNGK | ENSTIPDQRT | PVASTHSISS | AATPDRIKRP | 540 |
| 541 | RGTASRSTFH | GQPRERRTAT | YNGPPASPSL | SHEATPLSQT | RSRGSTNLF | KLTSKLTISR | 600 |
| 601 | NVSAEQKDEN | KEAKPRSLRF | TWSMKTSSM | DPGDMREIR | KVLDANNCY | EQRRERLLFC | 660 |
| 661 | VHGDGHAENL | VQWEMEVC | KL | PRLSLNGVRF | KRISGTSIAF | KNIASKIANE | 720 |

blue: MARK3 sequence expressed in fusion protein

¹NCBI/Protein accession number NP_001122392.2