

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

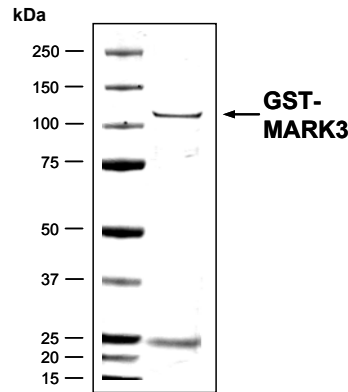
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

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

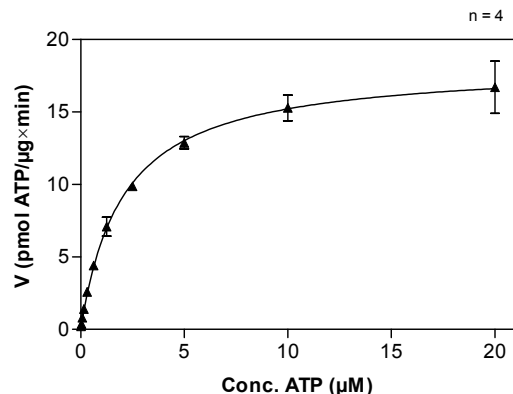
ATP-K_M: 2 µM

**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 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



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	DFLSKLPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHGG	RRRASVAAGI	240
241	LVPRGSPGLD	GIYARQMSTR	TPLPTVNERD	TENHTSHGDG	RQEVTSRTSR	SGARCRNSIA	300
301	SCADEQPHIG	NYRLLKTIGK	GNFAKVKLAR	HILTGREVAI	KIIDKTQLNP	TSLQKLFREV	360
361	RIMKILNHPN	IVKLFVIET	EKTLYLIMEY	ASGGEVFDYL	VAHGRMKEKE	ARSKFRQIVS	420
421	AVQYCHQKRI	VHRDLKAENL	LLDADMNIKI	ADFGFSNEFT	VGGKLDTFCG	SPPYAAPELF	480
481	QGKKYDGPEV	DVWSLGVILY	TLVSGSLPFD	GQNLKELRER	VLRGKYRIPF	YMSTDCENLL	540
541	KRFLVLNPIK	RGTLEQIMKD	RWINAGHEED	ELKPFVEPEL	DISDQKRIDI	MVGMGYSQEE	600
601	IQESLSKMKY	DEITATYLLL	GRKSSEVRPS	SDLNNSTGQS	PHHKVQRSVS	SSQKQRRYS	660
661	HAGPAIPSVV	AYPKRSQTST	ADSDLKEDGI	SSRKSSGS	SAVGGKIAPAS	MLGNASPNK	720
721	ADIPERKSS	TPSSNTASG	GMTRRNTYVC	SERTTADRHS	VIQNGKENST	IPDQRTPVAS	780
781	THSISSAATP	DRIRFPRGTA	SRSTFHGQPR	ERRTATYNGP	PASPPLSHEA	TPLSQTRSRG	840
841	STNLFSLKLS	KLTRSRNVSA	EQKDENEKAK	PRSLRFTWSM	KTTSSMDPGD	MMREIRKVL	900
901	ANNCYEQRE	RFLLCVHGD	GHAENLVQWE	MEVCKLPRLS	LNGVRFKRIS	GTSIAFKNIA	960
961	SKIANELKL						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	PHIGNYRLLK	60
61	TIGKGNFAKV	KLARHILTGR	EVAIKIIDKT	QLNPTSLQKL	FREVRIMKIL	NHPNIVKLFE	120
121	VIETEKTLYL	IMEYASGGEV	FDYLVAHGRM	KEKEARSKFR	QIVSAVQYCH	QKRIVHRDLK	180
181	AENLLLDADM	NIKIADFGFS	NEFTVGGKLD	TFCGSPPYAA	PELFQGKKYD	GPEVDVWSLG	240
241	VILYTLVSGS	LPFDGQNLKE	LRERVLRGKY	RIPFYMSTDC	ENLLKRFLVL	NPIKRGTLEQ	300
301	IMKDRWINAG	HEEDELKPFV	EPELDIDSDQK	RIDIMVGMGY	SQEEIQESLS	KMKYDEITAT	360
361	YLLGRKSSE	VRPSSDLNNS	TGQSPHHKVQ	RSVSSSQQR	RYSDHAGPAI	PSVVAYPKRS	420
421	QTSTADSDLK	EDGISSRKSS	GSAVGGKIA	PASPMLGNAS	NPNKADIPER	KKSSTVPSSN	480
481	TASGGMTRRN	TYVCSERTTA	DRHSVIQNGK	ENSTIPDQRT	PVASTHSISS	AATPDRIKRF	540
541	RGTASRSTFH	GQPRERTAT	YNGPPASPSL	SHEATPLSQT	RSRGSTNLF	KLTSKLTRSR	600
601	NVSAEQDEN	KEAKPRSLRF	TWSMKTSSM	DPGDMREIR	KVLDANNCYD	EQRRFLLFC	660
661	VHGDGHAENL	VQWEMEVCKL	PRLSLNGVRF	KRISGTSIAF	KNIASKIANE	LKL	720

blue: MARK3 sequence expressed in fusion protein

¹NCBI/Protein accession number NP_001122392.2