

MEKK3

mitogen-activated protein kinase kinase kinase 3

Recombinant Human Active Protein Kinase

HGNC Symbol: MAP3K3

Synonyms: MAPKKK3, MEK kinase3

Product No.: 0859-0000-1

Lot: 005

Description: Human MEKK3, full length, amino acids M₁-Y₆₇₂ (as in [NCBI/Protein](#) entry NP_002392.2), N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: MEKK3 Lot 005, was confirmed as MEKK3 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 99,304 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

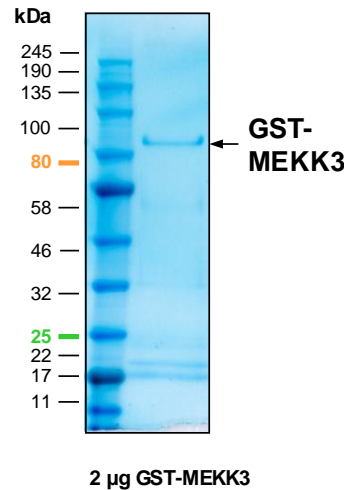
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.264 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

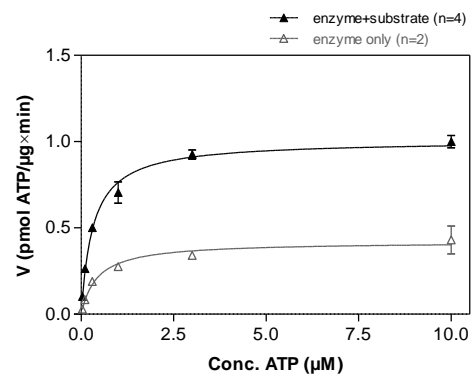
Biochemical Parameters:

Specific kinase activity (P_i transfer): 1 pmol/µg × min
ATP-K_M: 0.32 µM

MEKK3 Lot 005: Coomassie stain



MEKK3 Lot 005: 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, 20 µg/ml
 - Kinase: 4 µg/ml
- Filter binding assay
 - MSFC membrane (Millipore)

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.

MEKK3

Product No.: 0859-0000-1

GST-MEKK3 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVLFGQ	240
241	PLAMMDEQEA	LNSIMNDLVA	LQMNRRHRMP	GYETMKNKDT	GHSNRQSDVR	IKFEHNGERR	300
301	IIAFSRPVKY	EDVEHKVTV	FGQPLDLHYM	NNELSILLKN	QDDLK AIDI	LDRSSSMKSL	360
361	RILLLSQDRN	HNSSSPHSGV	SRQVRIKASQ	SAGDINTIYQ	PPEPRSRHLS	VSSQNPGRSS	420
421	PPPGYVPERQ	QHIAHQGSYT	SINSEGEFIP	ETSEQCMLDP	LSSAENSLSG	SCQSLDRSAD	480
481	SPSFRKSRMS	RAQSFDPNRQ	EYSDRETQLY	DKGVKGGTYP	RRYHVSVHHK	DYSDGRRTFP	540
541	RIRRHQGNLF	TLVPSSRSL	TNGENMGLAV	QYLDPRGRLR	SADSENALSV	QERNVPTKSP	600
601	SAPINWRRGK	LLQGQAFGRV	YLCYD VDTGR	ELASKQVQFD	PDSPETSKEV	SALECEIQLL	660
661	KNLQHERIVQ	YQGCLRDAE	KTLTIFMEYM	PGGSVKDQLK	AYGALTESVT	RKYTRQILEG	720
721	MSYLHSNMIV	HRDIKGANIL	RDSAGNVKLG	DFGASKRLQT	ICMSGTGMRS	VTGTPYWMSP	780
781	EVISGEGYGR	KADVWSLGCT	VVEMLTEKPP	WAEYEAMAAI	FKIATQPTNP	QLPSHISEHG	840
841	RDFLRRI FVE	ARQRPSAEEL	LTHHFAQLMY				900

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: MEKK3

MEKK3 wt ¹ Amino Acid Sequence							
1	MDEQEALNSI	MNDLVALQMN	RRHRMPGYET	MKNKDTGHSN	RQSDVRIKFE	HNGERRIAF	60
61	SRPVKYEDVE	HKVTTVFGQP	LDLHYMNNEL	SILLKNQDDL	DKAIDILDRS	SSMKSLRILL	120
121	LSQDRNHNS	SPHSGVSRQV	RIKASQSAGD	INTIYQPPPEP	RSRHLVSSQ	NPGRSSPPPG	180
181	YVPERQQHIA	RQGSYTSINS	EGEFIPETSE	QCMLDPLSSA	ENSLSGSCQS	LDRSADSPSF	240
241	RKSRMSRAQS	FDPNRQEYSD	RETQLYDKGV	KGGTYP RRYH	VSVHKKDYS	GRRTFPRIIR	300
301	HQGNLFTLVP	SSRSLSTNGE	NMGLAVQYLD	PRGRLRSADS	ENALSVQERN	VPTKSPSAPI	360
361	NWRRGKLLGQ	GAFGRVYLCY	DVDTGRELAS	KQVQFDPDSP	ETSKEVSALE	CEIQLLKNLQ	420
421	HERIVQYYGC	LRDRAEKTLT	IFMEYMPGGS	VKDQLKAYGA	LTESVTRKYT	RQILEGMSYL	480
481	HSNMIVHRDI	KGANILRDSA	GNVKLGDFGA	SKRLQ TICMS	GTGMRSVTGT	PYWMSPEVIS	540
541	GEGYGRKADV	WSLGCTV VEM	LTEKPPWAEY	EAMAAIFKIA	TQPTNPQLPS	HISEHGRDFL	600
601	RRIFVEARQR	PSAEELLTHH	FAQLMY				660

blue: MEKK3 sequence expressed in recombinant protein

¹NCBI/Protein accession number NP_002392.2