

ProQinase™ MKK3

mitogen-activated protein kinase kinase 3

Recombinant Human Active Protein Kinase

HGNC Symbol: MAP2K3

Synonyms: MEK3; MAPKK3; SAPKK2

Product No.: 1965-0000-1

Lot: 001

Description: Human MKK3, full length, amino acids M₁-S₃₄₇ (as in [NCBI/Protein](#) entry NP_659731.1), N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: MKK3 Lot 001, has been verified by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 69520 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: With MEKK3

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 10 µl are not advisable. Avoid repeated freeze-thaw cycles!

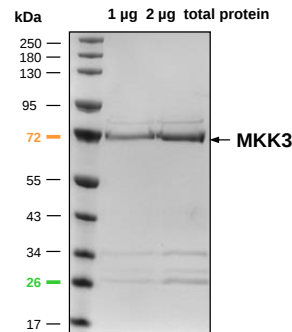
Protein concentration: 0.263 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

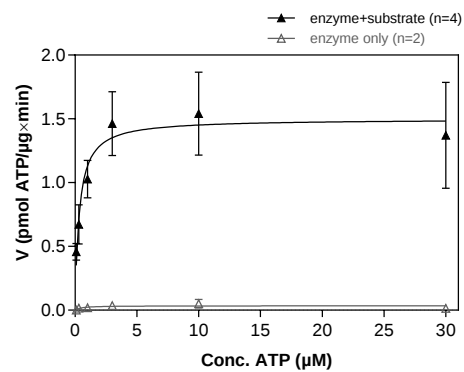
Specific kinase activity (P_i transfer): 1.5 pmol/µg*min

ATP-K_M: 0.33 µM

MKK3 Lot 001: Coomassie stain



MKK3 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: p38alpha K53A 40 µg/ml
 - Kinase: 4 µg/ml

Assay technology:
Radiometric filter binding assay
MSFC membrane (96 well plate, Millipore)

Sequence information

GST-MKK3 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLIERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPEML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHHG	RDSLEVLFGG	240
241	MESPASSQP	ASMPQSKGKS	KRKKDLRISC	MSKPPAPNPT	PPRNLDSTF	ITIGDRNFEV	300
301	EADDLVTISE	LGRGAYGVVE	KVRHAQSGTI	MAVKRIRATV	NSQEQKRLLM	DLINMRTVD	360
361	CFYTVTFYGA	LFREGDVWIC	MELMDTSLDK	FYRKVLDKNM	TIPEDILGEI	AVSIVRALEH	420
421	LHSKLSVIHR	DVKPSNVLIN	KEGHVKMCDF	GISGYLVDSV	AKTMDAGCKP	YMAPERINPE	480
481	LNQKGYNVKS	DVWSLGITMI	EMAILRFPYE	SWGTPFQQLK	QVVEEPSQPL	PADRFSPFEV	540
541	DFTAQCLRKN	PAERMSYLEL	MEHPFFTLHK	TKKTDIAAFV	KEILGEDSPW	ELAAACRVPP	600
601	GDLYRLC						660

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: MKK3

MKK3 wt ¹ Amino Acid Sequence							
1	MESPASSQPA	SMPQSKGKSK	RKKDLRISCM	SKPPAPNPTP	PRNLDSTFI	TIGDRNFEVE	60
61	ADDLVTISEL	GRGAYGVVEK	VRHAQSGTIM	AVKRIRATVN	SQEQKRLMD	LDINMRTVDC	120
121	FYTVTFYGA	FREGDVWICM	ELMDTSLDKF	YRKVLDKNMT	IPEDILGEIA	VSIVRALEHL	180
181	HSKLSVIHRD	VKPSNVLINK	EGHVKMCDFG	ISGYLVDSVA	KTMDAGCKPY	MAPERINPEL	240
241	NQKGYNVKSD	VWSLGITMIE	MAILRFPYES	WGTPFQQLKQ	VVEEPSQPLP	ADRFSPFEVD	300
301	FTAQCLRKNP	AERMSYLELM	EHPFFTLHKT	KKTDIAAFVK	EILGEDS		360

blue: MKK3 sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_659731.1