

MKK7

mitogen-activated protein kinase kinase 7

Recombinant Human Active Protein Kinase

HGNC Symbol: MAP2K7

Synonyms: MAPKK7, MEK7, PRKMK7

Product No.: 0726-0000-1

Lot: 004

Description: Human MKK7, full length, amino acids M₁-R₄₁₉ (as in [NCBI/Protein](#) entry NP_660186.1), N-terminal HIS₆ fusion protein with a Thrombin and TEV cleavage site, expressed in Sf9 insect cells

Product identity: MKK7 Lot 002, was confirmed as MKK7 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 52,809 Da Da

Expression host: Sf9 insect cells

Purification: Immobilized Metal Affinity Chromatography

Activation: With MEKK3

Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 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: 0.667 µg/µl

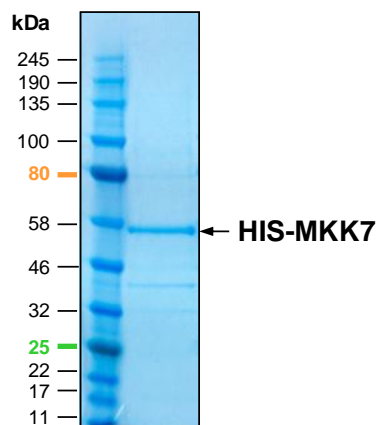
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

Specific kinase activity (P_i transfer): 12 pmol/µg × min

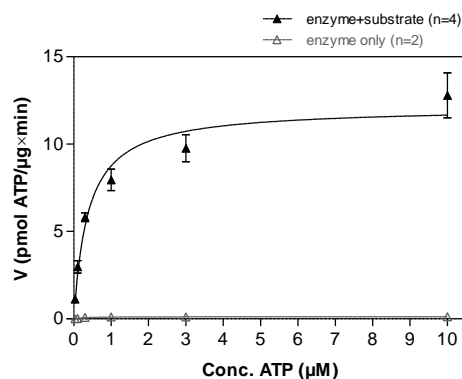
ATP-K_M: 0.4 µM

MKK7 Lot 004: Coomassie stain



1 µg HIS-MKK7 LOT004

MKK7 Lot 004: 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: JNK1 K55R/K56R 40 µg/ml
 - Kinase: 2 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.

MKK7

Product No.: 0726-0000-1

HIS-MKK7 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPIDPMG	HH	HHHGRRRAS	VAAGILVPRG	SPGLDGIYAR	TENLYFQGAM AASSLEQKLS	60
61	RLEAKLKQEN	REARRRIDLN	LDISPQRPRP	TLQLPLANDG	GSRSPSSESS	PQHPTPPARP	120
121	RHMLGLPSTL	FTPRSMESIE	IDQKLQEIMK	QTGYLTIGGQ	RYQAEINDLE	NLGEMGSGTC	180
181	GQVWKMFRK	TGHVIAVKQM	RRSGNKEENK	RILMDLDVVL	KSHDCPYIVQ	CFGTFITNTD	240
241	VFIAMELMGT	CAEKLKRMQ	GPIPERILGK	MTVAIVKALY	YLKEKHGVIH	RDVKPSNILL	300
301	DERGQIKLCD	FGISGRLVDS	KAKTRSAGCA	AYMAPERIDP	PDPTKPDYDI	RADVWSLGIS	360
361	LVELATGQFP	YKNCKTDFEV	LTKVLQEEPP	LLPGHMGFSG	DFQSFVKDCL	TKDHRKRPKY	420
421	NKLEHSFIK	RYETLEVDVA	SWFKDVMAKT	ESPRTSGVLS	QPHLPFFR		480

Red: HIS6-tag Pink: Thrombin cleavage site Green: TEV cleavage site blue: MKK7

MKK7 wt ¹ Amino Acid Sequence							
1	MAASSLEQKL	SRLEAKLKQE	NREARRRIDL	NLDISPQRPR	PTLQLPLAND	GGSRSPPSES	60
61	SPQHPTPPAR	PRHMLGLPST	LFTPRSMESI	EIDQKLQEIM	KQTGYLTIGG	QRYQAEINDL	120
121	ENLGEMSGT	CGQVWKMFRK	KTGHVIAVKQ	MRRSGNKEEN	KRILMDLDVV	LKSHDCPYIV	180
181	QCFGTFITNT	DVFIAMELMG	TCAEKLKRM	QGPIPERILG	KMTVAIVKAL	YYLKEKHGVI	240
241	HRDVKPSNIL	LDERGQIKLC	DFGISGRLVD	SKAKTRSAGC	AAYMAPERID	PPDPTKPDYD	300
301	IRADVWSLGI	SLVELATGQF	PYKNCKTDFE	VLTKVLQEEP	PLLPGHMGFS	GDFQSFVKDC	360
361	LTKDHRKRPK	YNKLEHSFI	KRYETLEVDV	ASWFKDVMAK	TESPRTSGVL	SQPHLPFFR	420

blue: MKK7 sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_660186.1

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.