

## ProQinase™ 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 M<sub>1</sub>-R<sub>419</sub> (as in [NCBI/Protein](#) entry NP\_660186.1), N-terminal HIS<sub>6</sub> 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**<sub>Fusion Protein</sub>: 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

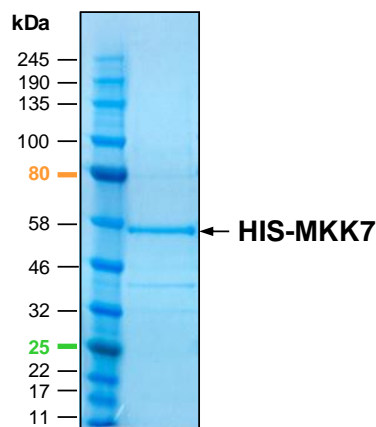
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Biochemical Parameters:**

Specific kinase activity (P<sub>i</sub> transfer): 12 pmol/µg × min

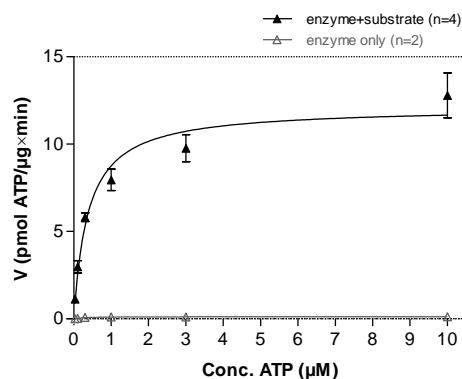
ATP-K<sub>M</sub>: 0.4 µM

### MKK7 Lot 004: Coomassie stain



1 µg HIS-MKK7 LOT004

### MKK7 Lot 004: Determination of V<sub>max</sub> and K<sub>M</sub> value for ATP



### Determination of K<sub>M</sub> value & Specific activity:

- Assay conditions:
  - 60 mM HEPES-NaOH, pH 7.5
  - 3 mM MgCl<sub>2</sub>
  - 3 mM MnCl<sub>2</sub>
  - 3 µM Na-orthovanadate
  - 1.2 mM DTT
  - 50 µg/ml PEG<sub>20,000</sub>
  - ATP (variable)
  - Substrate: JNK1 K55R/K56R 40 µg/ml
  - Kinase: 2 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

## ProQinase™ 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	GQVWKMRFRK	TGHVIAVKQM	RRSGNKEENK	RILMDLDVVL	KSHDCPYIVQ	CFGTFITNTD	240
241	VFIAMELMGT	CAEKLKRMQ	GPIPERILGK	MTVAIVKALY	YLKEKHGVIH	RDVKPSNILL	300
301	DERGQIKLCD	FGISGRIVDS	KAKTRSAGCA	AYMAPERIDP	PDPTKPDYDI	RADVWSLGIS	360
361	LVELATGQFP	YKNCKTDFEV	LTKVLQEEPP	LLPGHMGFSG	DFQSFVKDCL	TKDHRKRPKY	420
421	NKLLHSFIK	RYETLEVDVA	SWFKDVMAKT	ESPRTSGVLS	QPHLPFFR		480

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

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

blue: MKK7 sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](#) accession number NP\_660186.1