

## ProQinase™ MEK1 P124L

mitogen-activated protein kinase kinase 1

Recombinant Human Active Protein Kinase

HGNC Symbol: MAP2K1

Synonyms: MKK1, MAPKK1

Product No.: 1918-0000-1

Lot: 001

**Description:** Human MEK1, full length, amino acids M<sub>1</sub>-V<sub>393</sub> (as in [NCBI/Protein](#) entry NP\_002746.1), with a P124L point mutation, activated, N-terminal HIS-tag, expressed in Sf9 insect cells

**Product identity:** MEK1 P124L Lot 001, was confirmed as MEK1 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW**<sub>Fusion Protein</sub>: 46576 Da

**Expression host:** Sf9 insect cells

**Purification:** Immobilized Metal Affinity Chromatography

**Activation:** with RAF1

**Storage buffer:** 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 20 % glycerol

**Storage temperature:** -80°C

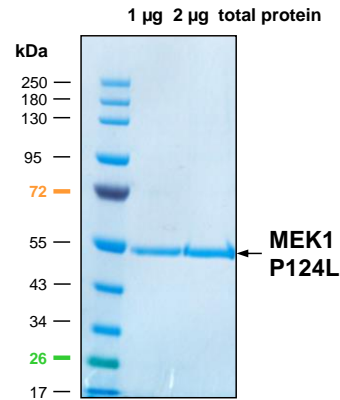
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

**Protein concentration:** 0.424 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

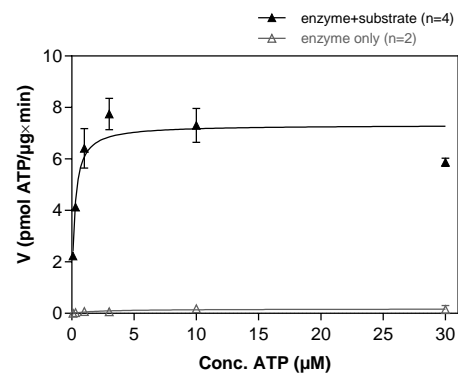
**Biochemical Parameters:**

Specific kinase activity (P<sub>i</sub> transfer): 7.3 pmol/µg x min  
ATP-K<sub>M</sub>: 0.2 µM

MEK1 P124L Lot 001:  
Coomassie stain



MEK1 P124L Lot 001:  
Determination of V<sub>max</sub> and K<sub>M</sub> value for ATP



- 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: ERK2 K54R 100 µg/ml
  - Kinase: 1 µg/ml
- Filter binding assay  
MSFC membrane (Millipore)

Recombinant Proteins

Sequence information

HIS-MEK1 P124L Recombinant Fusion Protein Amino Acid Sequence								
1	MSPILGDPMG	HHHHHH	GRDS	LEVLFGQ	PMP	KKKPTPIQLN	PAPDGSVAVNG TSSAETNLEA	60
61	LQKKLEEL	DEQQRK	RLEA	FLTQKQ	KVGE	LKDDDFEKIS	ELGAGNGGVV FKVSHKPSGL	120
121	VMARKLIHLE	IKPAIRNQII	RELQVLHECN	SIYIVGFYGA	FYSDGEISIC	MEHMDGGSLD		180
181	QVLKKAGRIP	EQILGKVSIA	VIKGLTYLRE	KHKIMHRDVK	PSNILVNSRG	EIKLCDFGVS		240
241	GQLIDSMANS	FVGTRSYMSP	ERLQGTHYSV	QSDIWSMGLS	LVEMAVGRYP	IPPPDAKELE		300
301	LMFGCQVEGD	AAETPPRPRT	PGRPLSSYGM	DSRPPMAIFE	LLDYIVNEPP	PKLPSGVFSL		360
361	EFQDFVNKCL	IKNPAERADL	KQLMVHAFIK	RSDAEEVDFA	GWLCSTIGLN	QPSTPTHAAG		420
421	V							480

Red: HIS6-tag Green: 3C cleavage site blue: MEK1 boxed: P124L mutation

MEK1 wt <sup>1</sup> Amino Acid Sequence								
1	MPKKKPTPIQ	LNPAPDGS	AV	NGTSSAETNL	EALQKKLEEL	ELDEQQRKRL	EAFLTQKQKV	60
61	GELKDDDFEK	ISELGAGNGG	VVFKVSHKPS	GLVMARKLIH	LEIKPAIRNQ	IIRELQVLHE		120
121	CNSYIVGFY	GAFYSDGEIS	ICMEHMDGGS	LDQVLKKAGR	IPEQILGKVS	IAVIKGLTYL		180
181	REKHKIMHRD	VKPSNILVNS	RGEIKLCDFG	VSQQLIDSMA	NSFVGTRSYM	SPERLQGTHY		240
241	SVQSDIWSMG	LSLVEMAVGR	YPIPPPD	AKE	LELMFGCQVE	GDAETPPRP	RTPGRPLSSY	300
301	GMDSRPPMAI	FELLDYIVNE	PPPKLPSGVF	SLEFQDFVNK	CLIKNPAERA	DLKQLMVHAF		360
361	IKRSDAEEVD	FAGWLCSTIG	LNQPSTP	THA	AGV			420

blue: kinase sequence expressed in recombinant protein Red: variant in recombinant protein

<sup>1</sup>NCBI/Protein accession number NP\_002746.1