

## ProQinase™ MEK1 K97M

mitogen-activated protein kinase kinase 1

### Recombinant Protein Kinase Substrate

HGNC Symbol: MAP2K1

Synonyms: PRKMK1, MAPKK1, MKK1

Product No.: 0785-0000-1

Lot: 075

**Description:** Human MEK1, full length, amino acids M<sub>1</sub>-V<sub>393</sub> (as in [NCBI/Protein](#) entry NP\_002746.1), mutationally inactivated by amino acid exchange K<sub>97</sub>M, untagged, expressed in Sf9 insect cells

**Theoretical MW**<sub>Fusion Protein</sub>: 43,784 Da

**Expression host:** Sf9 insect cells

**Purification:** GST-Affinity Chromatography

**ATPase activity:** In an ADP-Glo™ assay (Promega), the ATP → ADP conversion within 30 min is approx. 1% at a concentration of 100 µg/ml substrate.

Detailed ATPase assay conditions on request

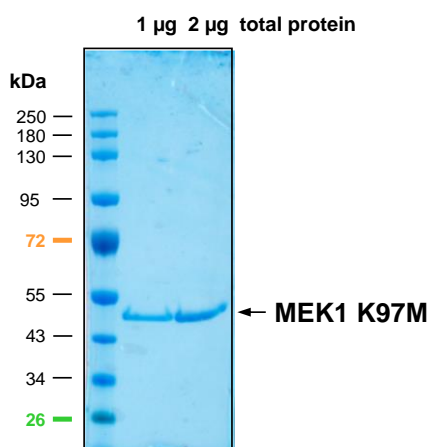
**Storage buffer:** 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 10 % glycerol

**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!

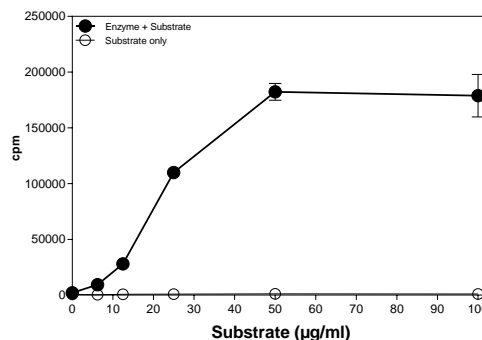
**Protein concentration:** 0.461 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**MEK1 K97M Lot 075:**  
**Coomassie stain**



### Phosphorylation of MEK1 K97M by RAF1 YDYD

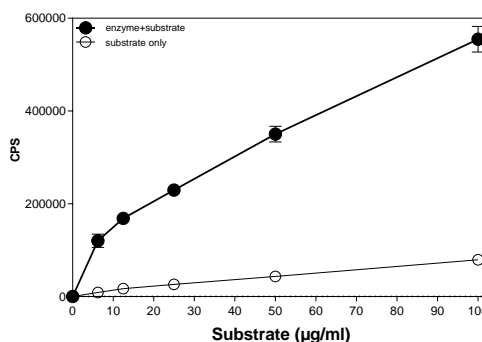
#### Radiometric filter binding assay



#### Assay conditions:

70 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: 1 µM  
Substrate: variable concentration  
Kinase: 1 µg/ml  
MSFC filter plates (Corning)

### ADP-Glo™ assay (Promega)



#### Assay conditions:

70 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: 10 µM  
1 % (v/v) DMSO  
Substrate: variable concentration  
Kinase: 2 µg/ml

MEK1 K97M

Product No.: 0785-0000-1

MEK1 K97M Recombinant Fusion Protein Amino Acid Sequence							
1	GPLAMTKKKP	TPIQLNPAPD	GSVNGTSSA	ETNLEALQKK	LEEELEDEQQ	RKRLEAFLTQ	60
61	KQKVGELKDD	DFEKISELGA	GNGGVVFKVS	HKPSGLVMAR	MLIHLEIKPA	IRNQIIRELQ	120
121	VLHECNSPYI	VGFGAFYSD	GEISICMEHM	DGGSLDQVLK	KAGRIPEQIL	GKVSIAVIKG	180
181	LTYLREKHKI	MHRDVKPSNI	LVNSRGEIKL	CDFGVSGQLI	DSMANSFVGT	RSYMSPERLQ	240
241	GTHYSVQSDI	WSMGLSLVEM	AVGRYPIPPP	DAKELELMFG	CQVEGDAAET	PPRPRTPGRP	300
301	LSSYGMDSRP	PMAIFELLDY	IVNEPPPKLP	SGVFSLEFQD	FVNKCLIKNP	AERADLKQLM	360
361	VHAFIKRSDA	EEVDFAGWLC	STIGLNQPST	PTHAAGV			420

1-4: legacy of tag cleavage    blue: MEK1    boxed: variation from RefSeq

This product was manufactured at Reaction Biology in Freiburg, Germany, and is for in vitro research use only, not for use in human or animals.  
© European Union, 2020. Material may not be reproduced or distributed without written permission from Reaction Biology Europe GmbH