

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: 070

Description: Human MEK1, full length, amino acids M₁-V₃₉₃ (as in [NCBI/Protein](#) entry NP_002746.1), mutationally inactivated by amino acid exchange K₉₇M, untagged, expressed in Sf9 insect cells/E.coli

Theoretical MW_{Fusion Protein}: 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

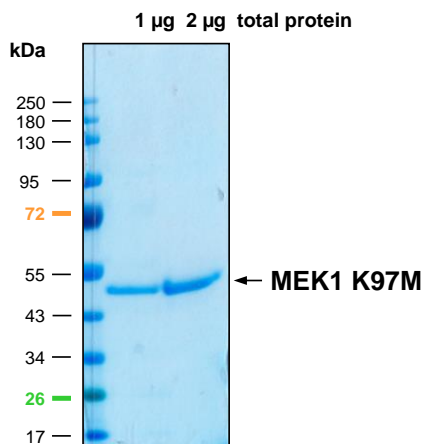
Storage temperature: -80°C

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

Protein concentration: 0.464 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

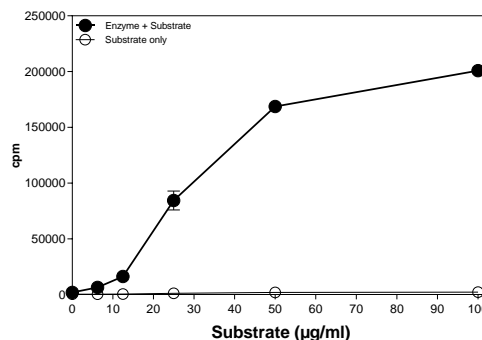
Substrate Lot 070:

Coomassie stain



Phosphorylation of SUBSTRATE by KINASE

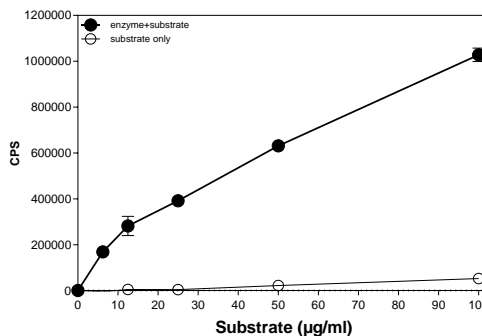
Radiometric filter binding assay



Assay conditions:

70 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: 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 MgCl₂
3 mM MnCl₂
3 µM Na-orthovanadate
1.2 mM DTT
50 µg/ml PEG_{20,000}
ATP: variable concentration
1 % (v/v) DMSO
Substrate: 100 µg/ml
Kinase: 2 µg/ml

MEK1 K97M

Product No.: 0785-0000-1

| MEK1 K97M Recombinant Fusion Protein Amino Acid Sequence | | | | | | | |
|--|------------|------------|------------|------------|------------|------------|-----|
| 1 | GPLAMTKKKP | TPIQLNPAPD | GSAVNGTSSA | ETNLEALQKK | LEELELDEQQ | RKRLEAFLTQ | 60 |
| 61 | KQKVGELKDD | DFEKISELGA | GNGGVVFKVS | HKPSGLVMAR | MLIHLEIKPA | IRNQIIRELQ | 120 |
| 121 | VLHECNSPYI | VGFYGFYSD | GEISICMEHM | DGGSLDQVLK | KAGRIPEQIL | GKVSIAVIKQ | 180 |
| 181 | LTYLREKHKI | MHRDVKPSNI | LVNSRGEIKL | CDFGVSGQLI | DSMANSFVGT | RSYMSPERLQ | 240 |
| 241 | GTHYSVQSDI | WSMGLSLVEM | AVGRYPIPPP | DAKELELMFG | CQVEGDAEET | PPRPRTGGRP | 300 |
| 301 | LSSYGMDSRP | PMAIFELLDY | IVNEPPPPLP | 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