

Certificate of Analysis

MAPKAP Kinase 2, active

(Recombinant enzyme expressed in *E.coli* cells)

Item # 14-337, 14-337-K, 14-337M

Parent Lot # WAE0265

The data presented in this document apply to the parent lot shown above and to all pack sizes derived from subsequent vialling runs of this parent lot. An alphabetical suffix after the parent lot number is used to denote each vialling run.

Product Description: N-Terminal GST-tagged, recombinant, human MAPKAP Kinase 2, amino acids 46–end, expressed in *E.coli* cells. Purified using glutathione agarose. Activated using MAP Kinase 2 and repurified using Q-sepharose. Purity 84% by SDS-PAGE and Coomassie blue staining. MW = 70.2kDa.

Formulation: 0.77mg/ml of enzyme in 50mM Tris/HCl, 150mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 50% glycerol, 1mM benzamidine, 0.2mM PMSF, 0.1% 2-mercaptoethanol, 5mM dithiothreitol. Liquid at -20°C.

Specific Activity (Parent lot# WAE0265): 3002U/mg, where one unit of MAPKAP Kinase 2, active activity is defined as 1nmol phosphate incorporated into 30µM substrate peptide (KKLNRTLVA, cat# 12-240) per minute at 30°C with a final ATP concentration of 100µM.

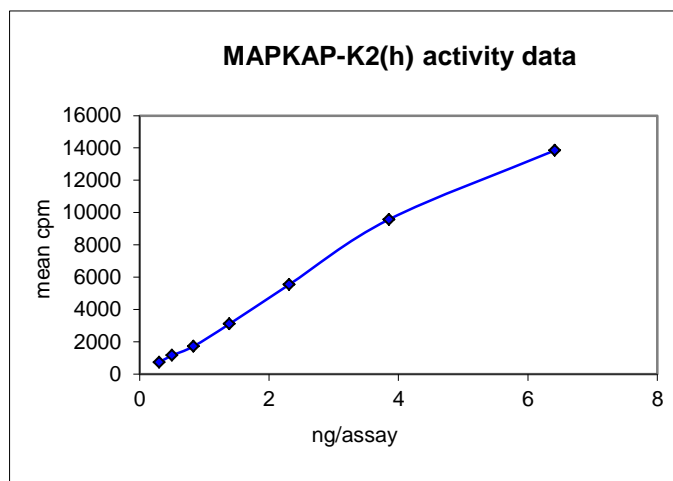
Storage and Stability: On receipt of material store at -20°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

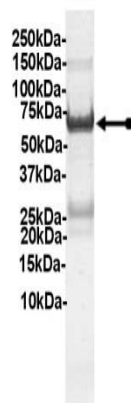
Quality Control Testing

Kinase Assay: 0.30–6.42ng of this lot of enzyme phosphorylated 30µM substrate peptide (KKLNRTLVA) in the assay described on page two. Assay background was subtracted from the actual counts to yield the results shown below.

MS Tryptic Fingerprint: Confirmed product identity as MAPKAP Kinase 2 with the translated native sequence listed on page three.



SDS-PAGE and Coomassie Stain: Purity was assessed by SDS-PAGE and Coomassie blue staining using 3µg of MAPKAP Kinase 2, active.



Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

- 1. 20 x Reaction Buffer:** 1M Na- β -glycerol phosphate, 2mM EGTA pH7.5.
- 2. MAPKAP Kinase 2 substrate peptide (KKLNRTLVA Catalogue# 12-240):** Use at a final assay concentration of 30 μ M. Make a 300 μ M stock. Add 2.5 μ l of stock per assay point.
- 3. MAPKAP Kinase 2, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 5% glycerol, 0.01% Brij-35, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 0.30–6.42ng per assay point.
- 4. [γ -³³P]ATP:** 2.5 x magnesium acetate/[γ -³³P]ATP cocktail: 25mM MgAc and 0.25mM ATP to which is added [γ -³³P]ATP (specific activity approximately 500 - 800cpm/pmol as required.)

Assay Procedure (96 well plate format):

1. Add 5 μ l of 20 x reaction buffer per well.
2. Add 2.5 μ l of substrate peptide (**KKLNRTLVA**).
3. Add **2.5 μ l (0.30–6.42ng) MAPKAP-K2, active**.
4. Add 5 μ l of dH₂O.
5. Add 10 μ l of diluted [γ -³³P]ATP mixture.
6. Incubate for 10 minutes at 30°C.
7. Stop the reaction by adding 5 μ l of 3% phosphoric acid.
8. Transfer a 10 μ l aliquot onto the appropriate area of a **P30 Filtermat**.
9. Wash the filtermat three times for 5 minutes with 50mM phosphoric acid.
10. Wash the filtermat once for 2 minutes with methanol.
11. Transfer the filtermat to a sealable plastic bag and add 4ml of scintillation cocktail.
12. Read in a scintillation counter. Compare cpm of enzyme samples with cpm of control samples that contain all assay components plus 1 μ l of 30% phosphoric acid.

Certificate of Analysis

MAPKAP Kinase 2 Sequence Information

<u>Protein</u>	human MAPKAP Kinase 2
<u>Tags</u>	N-Terminal GST
<u>Native sequence</u>	F243 of the recombinant protein is equivalent to F46 of human MAPKAP Kinase 2
<u>Accession number</u>	GenBank NM_032960. The recombinant sequence is D116H (native coordinates) with respect to GenBank NM_032960. This conflict is reported in SWISSPROT P49137. The cDNA also contains the translational conflict A399G, and a c-myc epitope at the C-terminus.

Recombinant MAPKAP Kinase 2 amino acid sequence:

```

1  MSPILGYWKI  KGLVQPTRL  LEYLEEKYEE  HLYERDEGDK  WRNKKFELGL  EFPNLPYYID
61  GDVKLTQSMA  IIRYIADKHN  MLGGCPKERA  EISMLEGAVL  DIRYGVSRIA  YSKDFETLKV
121  DFLSKLP EML  KMFEDRLCHK  TYLNGDHVTH  PDFMLYDALD  VVLYMDPMCL  DAFPKLVCFK
181  KRIEAI PQID  KYLKSSKYIA  WPLQGWQATF  GGGDHPPKSD  LVPRGSPGIS  GGGGGILEAT
241  MEFHVK SGLQ  IKKNAIDDDY  KVTSQVLGLG  INGKVLQIFN  KRTQEKFALK  MLQDCPKARR
301  EVELHW RASQ  CPHIVRIVDV  YENLYAGRKC  LLIVMECLDG  GELFSRIQDR  GDQAFTEREA
361  SEIMKSI GEA  IQYLHSINIA  HRDVKPENLL  YTSKRPNAIL  KLTDFGFAKE  TTSHNSLTTP
421  CYTPYYVA PE  VLGPEKYDKS  CDMWSLGVIM  YILLCGYPPF  YSNHGLAISP  GMKTRIRMGQ
481  YEFNP EWSE  VSEEVKMLIR  NLLKTEPTQR  MTITEFMNHP  WIMQSTKVPQ  TPLHTRSRLK
541  EDKERW EDVK  EEMTSALATM  RVDYEQIKIK  KIEDASNPLL  LKRRKKARAL  EAAALGHMEQ
601  KLISEEDLK

```

Recombinant MAPKAP Kinase 2 nucleotide sequence:

```

1  atgtccccta tactaggtta ttggaaaatt aagggccttg tgcaaccac tcgacttctt
61  ttggaatata ttgaagaaaa atatgaagag catttgtatg agcgcgatga aggtgataaa
121  tggcgaagaa aaaagtttga attgggtttg gagtttccca atcttcctta ttatattgat
181  ggtgatgtta aattaacaca gtctatggcc atcatagctt atatactgta caagcacaac
241  atgttgggtg gttgtccaaa agagcgtgca gagatttcaa tgcttgaagg agcggttttg
301  gatattagat acggtgtttc gagaattgca tatagtaaag actttgaaac tctcaaagtt
361  gattttctta gcaagctacc tgaatgctg aaaatgttcg aagatcgttt atgtcataaa
421  acatatttaa atggtgatca tgtaaccat cctgacttca tgttgtatga cgctcttgat
481  gttgttttat acatggacc aatgtgcctg gatgcgttcc caaaattagt ttgttttaaa
541  aaacgtattg aagctatccc acaaattgat aagtacttga aatccagcaa gtatatagca
601  tggcctttgc agggctggca agccacgttt ggtggtggcg accatcctcc aaaatcggat
661  ctggttccgc gtggatcccc ggaatttcc ggtggtggcg gtggaattct agaggccacc
721  atggagtacc acgtcaagtc cggcctgcag atcaagaaga acgcatcatc cgatgactac
781  aaggtcacca gccaggtcct ggggctgggc atcaacggca aagttttgca gatcttcaac
841  aagaggacc caggagaaat cgccctcaaa atgcttcagg actgccccaa ggcccgcagg
901  gaggtggagc tgactggcg ggcctcccag tgcccgcaca tcgtacggat cgtggatgtg
961  tacgagaatc tgtacgcagg gaggaagtgc ctgctgattg tcatggaatg tttggacggt
1021 ggagaactct ttagcgaat ccaggatcga ggagaccagg cattcacaga aagagaagca
1081 tccgaaatca tgaagagcat cggtagggcc atccagtatc tgcattcaat caacattgcc
1141 catcgggatg tcaagcctga gaatctctta tacacctcca aaaggcccaa cgccatcctg
1201 aaactcactg actttggctt tgccaaggaa accaccagcc acaactctt gaccactcct
1261 tgttatacac cgtactatgt ggctccagaa gtgctgggtc cagagaagta tgacaagtcc
1321 tgtgacatgt ggtccctggg tgtcatcatg tacatcctgc tgtgtgggta tcccccttc
1381 tactccaacc acggccttgc catctctccg ggcattgaaga ctgcacatcc aatgggcccag

```

Certificate of Analysis

```
1441 tatgaatttc ccaaccaga atggtcagaa gtatcagagg aagtgaagat gtcattcgg
1501 aatctgctga aaacagagcc caccagaga atgaccatca ccgagtttat gaaccaccct
1561 tggatcatgc aatcaacaaa ggtccctcaa acccactgc acaccagccg ggtcctgaag
1621 gaggacaagg agcgggtgga ggatgtcaag gaggagatga ccagtgcctt ggccacaatg
1681 cgcgttgact acgagcagat caagataaaa aagattgaag atgcatccaa ccctctgctg
1741 ctgaagaggc ggaagaaagc tcgggccttg gaggctgcgg ctctgggcca catggagcag
1801 aagctgatca gcgaggagga cctgaagtga
```

Reviewed and approved by site quality representative.

Unless otherwise stated in our catalogue or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

© 2014 **Eurofins Pharma Discovery Services UK Limited** is an independent member of Eurofins Discovery Services