

## Certificate of Analysis

### B-Raf (V599E), active

#### Also designated B-Raf (V600E)

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-557, 14-557-K, 14-557M

Parent Lot # 28439U

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 B-Raf residues 416–end, containing a V599E mutation. Expressed by baculovirus in Sf21 insect cells. Purified using glutathione agarose. Purity 44% by SDS-PAGE and Coomassie blue staining. MW = 67.3kDa.

**Specific Activity (Parent lot# 28439U):** 167914U/mg, where one unit of B-Raf, active = 1 unit of MAPK2(m) activity which in turn is equivalent to 1nmol phosphate incorporated into 0.33mg/ml myelin basic protein per minute at 30°C with a final ATP concentration of 100µM. Note the activity is determined by a triple linked assay which involves the activation of MEK1(h) by B-Raf, followed by the subsequent activation of MAPK2(m) by the activated MEK1(h).

**Formulation:** 0.823mg/ml of enzyme in 50mM Tris/HCl pH7.5, 150mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 1mM benzamidine, 0.2mM PMSF, 0.1% 2-mercaptoethanol. Frozen solution.

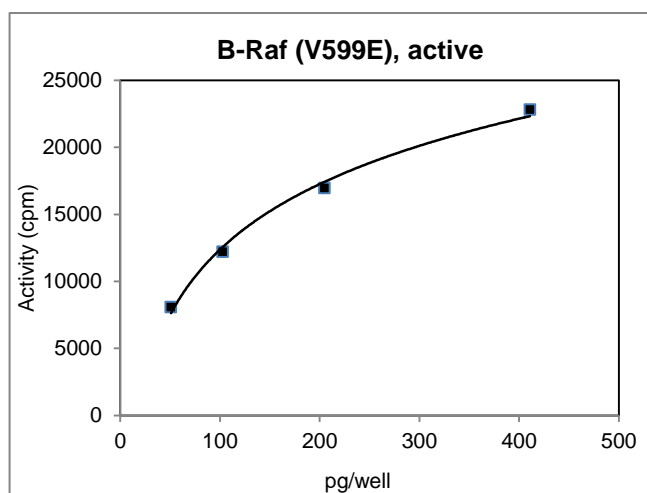
**Storage and Stability:** On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

**Handling Recommendations:** Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled micro-centrifuge tubes and immediately snap-freeze the vials in liquid nitrogen prior to re-storage at -70°C.

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

### Quality Control Testing

**Kinase Assay:** 51-411pg of this lot of enzyme was used to activate 0.2µM MEK1(h) in the assay described on page two. Assay background was subtracted from the actual counts to yield the results.



**MS Tryptic Fingerprint:** Confirmed product identity as B-Raf with the translated sequence listed on page three.



## Certificate of Analysis

### Kinase Assay Protocol

#### Stock Solutions:

- 1. 10 x Reaction Buffer:** 500mM Tris/HCl pH7.5, 1mM EGTA, 1mM Na<sub>3</sub>VO<sub>4</sub>, 1% 2-mercaptoethanol, 0.1% Brij-35.
- 2. MEK1(h), unactive:** Use at a final assay concentration of 0.2μM (0.0126mg/ml). Prepare a 0.126mg/ml stock and add 2.5μl stock per assay point.
- 3. MAPK2(m), unactive:** Use at a final assay concentration of 2μM (0.14mg/ml). Prepare a 0.7mg/ml stock and add 5μl of stock per assay point.
- 4. Myelin Basic Protein (MBP):** Use at a final assay concentration of 0.33mg/ml. Prepare a 3.33mg/ml stock and add 2.5μl stock per assay point.
- 5. B-Raf, active:** Dilute with 25mM Tris/HCl pH7.5, 0.1mM EGTA, 1mg/ml BSA. Use 51-411pg per assay point.
- 6. [γ-<sup>33</sup>P]ATP:** 2.5 x magnesium acetate/[γ-<sup>33</sup>P]ATP cocktail: 25mM MgAc and 0.25mM ATP to which is added [γ-<sup>33</sup>P]ATP (specific activity approximately 500 - 800cpm/pmol as required.)

#### Assay Procedure (96 well plate format):

1. Add 2.5μl of 10 x reaction buffer to wells.
2. Add 2.5μl of **MEK1(h), unactive**.
3. Add 2.5μl of **MAPK2 (m), unactive**.
4. Add 2.5μl of **MBP**.
5. Add **5μl (51–411pg) B-Raf, active**.
6. Add 10μl of diluted [γ-<sup>33</sup>P] ATP mixture.
7. Incubate for 10 minutes at 30°C.
8. Stop the reaction by adding 5μl 3% phosphoric acid to each well.
9. Using a multichannel pipette, spot 10μl onto the appropriate area of a **P30 Filtermat**.
10. Wash the filtermat twice for 5 minutes with 75mM phosphoric acid.
11. Wash the filtermat once for 2 minutes with methanol.
12. Once dry, transfer the filtermat to a sealable plastic bag and add 4ml of scintillation cocktail.
13. Read in a scintillation counter. Compare cpm of enzyme samples with cpm of control samples that contain all assay components plus 1μl of 30% phosphoric acid.

## Certificate of Analysis

### B-Raf (V599E) Sequence Information

<b><u>Protein</u></b>	Human B-Raf (V599E)
<b><u>Tags</u></b>	N-Terminal GST
<b><u>Native sequence</u></b>	Q237 of the recombinant protein is equivalent to Q416 of human B-Raf
<b><u>Accession number</u></b>	GenBank NM_004333. The V599E mutation is thought to mimic phosphorylation of the native enzyme, resulting in a protein with high activity and leading to constitutive ERK activation. As B-Raf is commonly activated by somatic point mutation in human cancer, it may provide a new therapeutic approach to malignant melanoma. (Davies H., <i>et al.</i> , <i>Nature</i> , 2002. <b>417</b> : 949-54, and Mercer KE., & Pritchard CA., <i>Biochim Biophys Acta</i> . 2003. <b>1653</b> : 25-40)

#### Recombinant B-RAF (V599E) amino acid sequence:

```

1  MSPILGYWKI  KGLVQPTRLL  LEYLEEKYEE  HLYERDEGDK  WRNKKFELGL  EFPNLPYYID
61  GDVKLTQ SMA  IIRYIADKHN  MLGGCPKERA  EISMLEGAVL  DIRYGVSRIA  YSKDFETLKV
121  DFLSKLP EML  KMFEDRLCHK  TYLNGDHVTH  PDFMLYDALD  VVLYMDPMCL  DAFPKLVCFK
181  KRIEAI P QID  KYLKSSKYIA  WPLQG WQATF  GGGDHPPKSD  LEVLFQGP EF  KGLRRQ QKSP
241  GPQRE R KSSS  SSEDNRN MKT  LGRRDSSDDW  EIPDGQITVG  QRIGSGSFGT  VYKKGWHGDV
301  AVKMLN V TAP  TPQQLQAFKN  EVGVL R KTRH  VNILLFMGYS  TKPQLAIVTQ  WCEGSSLYHH
361  LHI I ETKFEM  IKLIDIARQT  A QGM DY L HAK  S I IHRDLKSN  NIFLHEDLTV  KIGDFGLATE
421  KSRWSG SHQF  EQLSGSILWM  APEVIRMQDK  NPYSFQSDVY  AFGIVLYELM  TGQLPYSNIN
481  NRDQII F MVG  RGYLSPDL SK  VRSNCPKAMK  RLMAECLKKK  RDERPLFPQI  LASIELLARS
541  LPKIHR SASE  PSLNRAGFQT  EDFSLYACAS  PKTPIQAGGY  GAFPVH

```

#### Recombinant B-RAF (V599E) nucleotide sequence:

```

1  atgtccccta  tactaggtta  ttggaaaatt  aagggccttg  tgcaaccac  tcgacttctt
61  ttggaatata  ttgaagaaaa  atatgaagag  catttgtatg  agcgcgatga  aggtgataaa
121  tggcgaaaaca  aaaagtttga  attgggtttg  gagtttccca  atcttcctta  ttatattgat
181  ggtgatgtta  aattaacaca  gtctatggcc  atcatacggt  atatagctga  caagcacaac
241  atgttgggtg  gttgtccaaa  agagcgtgca  gagatttcaa  tgcttgaagg  agcggttttg
301  gatattagat  acggtgtttc  gagaattgca  tataagtaaag  actttgaaac  tctcaaagtt
361  gattttctta  gcaagctacc  tgaaatgctg  aaaatgttcg  aagatcgttt  atgtcataaa
421  acatatttaa  atggtgatca  tgtaaccat  cctgacttca  tgttgtatga  cgctcttgat
481  gttgttttat  acatggacc  aatgtgacct  gatgcgttcc  caaaattagt  ttgttttaa
541  aaacgtattg  aagctatccc  acaaattgat  aagtacttga  aatccagcaa  gtatatagca
601  tggcctttgc  agggctggca  agccacgttt  ggtggtggcg  accatcctcc  aaaatcggat
661  ctggaagtcc  tgttccaggg  gcccgaaatt  aaaggcctac  gtcgacaaca  gaaatctcca
721  ggacctcagc  gtgaaaggaa  gtcacttcca  tcctcagaag  acaggaatcg  aatgaaaaca
781  cttggtagac  gggactcgag  tgatgattgg  gagattcctg  atgggcagat  tacagtggga
841  caaagaattg  gatctggatc  atttggaaca  gtctacaagg  gaaagtggca  tgggtgatgtg
901  gcagtgaaaa  tgttgaatgt  gacagcacct  acacctcagc  agttacaagc  cttcaaaaat
961  gaagtaggag  tactcaggaa  aacacgacat  gtgaatatcc  tactcttcat  gggctattcc
1021  acaaagccac  aactggctat  tgttaccag  tgggtgtgagg  gctccagctt  gtatcaccat
1081  ctccatatca  ttgagaccaa  atttgagatg  atcaaactta  tagatattgc  acgacagact
1141  gcacagggca  tggattactt  acacgccaag  tcaatcatcc  acagagacct  caagagtaat
1201  aatatatttc  ttcatagaaga  cctcacagta  aaaataggtg  attttggctt  agctacagag
1261  aatctcagat  ggagtgggtc  ccatcagttt  gaacagttgt  ctggatccat  tttgtggatg
1321  gcaccagaag  tcatcagaat  gcaagataaa  aatccataca  gctttcagtc  agatgtatat
1381  gcatttggaa  ttgttctgta  tgaattgatg  actggacagt  taccttattc  aaacatcaac
1441  aacagggacc  agataatttt  tatggtggga  cgaggatacc  tgtctccaga  tctcagtaag
1501  gtacggagta  actgtccaaa  agccatgaag  agattaatgg  cagagtgcct  caaaaagaaa

```

## Certificate of Analysis

```
1561 agagatgaga gaccactctt tccccaaatt ctgcctcta ttgagctgct ggcccgctca
1621 ttgccccaaa ttcaccgcag tgcacagaa ccctccttga atcgggctgg tttccaaaca
1681 gaggatttta gtctatatgc ttgtgcttct ccaaaaacac ccatccaggc agggggatat
1741 ggtgcgtttc ctgtccactg a
```

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.