

## Certificate of Analysis

### FGFR1 (V561M), active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-734, 14-734-K, 14-734M

Parent Lot # D7BN045U

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 FGFR1 amino acids 456–765 containing the V561 mutation. Expressed by baculovirus in Sf21 insect cells. Purified using glutathione agarose. Co-ordinates encompass the protein kinase domain (478–767) according to SP P11362. The coordinates for recombinant human FGFR1 were described in Homann *et al*, J Biotech (2001); **86**:51-58. Blencke *et al* (2004) suggested mutations to a conserved threonine residue at the ATP binding site would result in inhibitor resistance. The amino acid valine 561 was mutated to a methionine in FGFR1 which corresponded to previously reported mutations found in Abl (T315) and EGFR (T766) which had been shown to confer resistance to selective inhibitors. Assay data for FGFR1 V561M had shown that this mutation had conferred resistance to PP58 (pyrido [2,3-d] pyrimidine tyrosine kinase inhibitor) compared with that of the wild type. (Blencke, S. *et al*, Chemistry & Biology (2004);**11**:691-701) Purity 100% by SDS-PAGE and Coomassie blue staining. MW = 62.3kDa.

**Specific Activity (Parent lot# D7BN045U):** 4741U/mg, where one unit of FGFR1 (V561M) activity is defined as 1nmol phosphate incorporated into 500 $\mu$ M (GGEEEEYFELVKKKK) per minute at 30°C with a final ATP concentration of 100 $\mu$ M.

**Formulation:** 1.2mg/ml of enzyme in 50mM Tris/HCl pH7.5, 300mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 1mM benzamidine, 0.2mM PMSF, 0.1% 2-mercaptoethanol. Frozen solution.

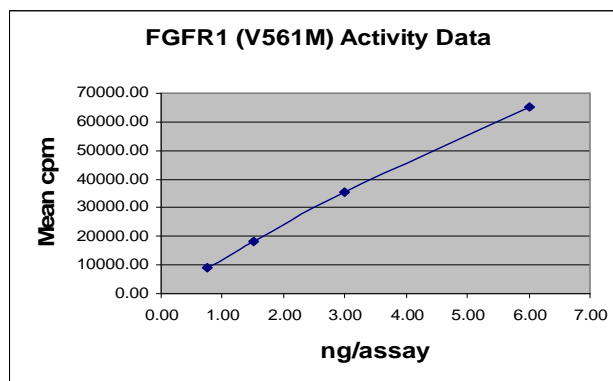
**Storage and Stability:** On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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

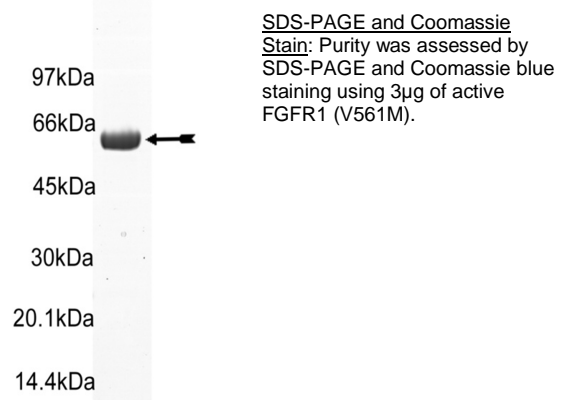
**FOR IN VITRO RESEARCH USE ONLY  
NOT FOR USE IN HUMANS OR ANIMALS**

### Quality Control Testing

**Kinase Assay:** 0.75–6.00ng of this lot of enzyme phosphorylated 500 $\mu$ M (GGEEEEYFELVKKKK) 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 FGFR1 (V561M) with the translated native sequence listed on page three.



## Certificate of Analysis

### Kinase Assay Protocol

#### Stock Solutions:

1. **5 x Reaction Buffer:** 40mM MOPS/NaOH pH7.0, 1mM EDTA.
2. **(GGEEEEYFELVKKKK):** Use at a final assay concentration of 500 $\mu$ M. Prepare a 5mM stock and add 2.5 $\mu$ l of stock per assay point.
3. **FGFR1 (V561M), active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 0.75–6.00ng per assay point.
4. **[ $\gamma$ -<sup>33</sup>P]ATP:** 2.5 x magnesium acetate/[ $\gamma$ -<sup>33</sup>P]ATP cocktail: 25mM MgAc and 0.25mM ATP to which is added [ $\gamma$ -<sup>33</sup>P]ATP (specific activity approximately 500 - 800cpm/pmol as required.)

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

1. Add 5 $\mu$ l of 5 x reaction buffer per assay to wells.
2. Add 2.5 $\mu$ l of **(GGEEEEYFELVKKKK)**.
3. Add **2.5 $\mu$ l (0.75–6.00ng) FGFR1 (V561M), active**.
4. Add 10 $\mu$ l of diluted [ $\gamma$ -<sup>33</sup>P]ATP mixture.
5. Add 5 $\mu$ l of dH<sub>2</sub>O.
6. Incubate for 10 minutes at 30°C.
7. Stop the reaction by adding 5 $\mu$ l of 3% phosphoric acid.
8. Transfer a 10 $\mu$ l aliquot onto the appropriate area of a **P30 Filtermat**.
9. Wash the filtermat three times for 5 minutes with 75mM 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 $\mu$ l of 30% phosphoric acid.

## Certificate of Analysis

### FGFR1 (V561M) Sequence Information

<b>Protein</b>	Human FGFR1 (V561M)
<b>Tags</b>	N-terminal GST
<b>Native sequence</b>	M231 of recombinant sequence is equivalent to M456 of native human FGFR1
<b>Accession number</b>	GenBank NM_000604

#### Recombinant FGFR1 (V561M) amino acid sequence:

```

1  MSPILGYWKI  KGLVQPTRL  LEYLEEKYEE  HLYERDEGDK  WRNKKFELGL  EFPNLPYYID
61  GDVKLTQSM  IIRYIADKHN  MLGGCPKERA  EISMLEGAVL  DIRYGVSRIA  YSKDFETLKV
121  DFLSKLP  EML  KMFEDRLCHK  TYLNGDHVTH  PDFMLYDALD  VVLYMDPMCL  DAFPKLVCFK
181  KRIEAI  PQID  KYLKSSKYIA  WPLQGWQATF  GGDHPPKSD  LEVLFQGF  EPEF  MLAGVSEYEL
241  PEDPRWEL  PR  DRLVLGKPLG  EGCFGQVVLA  EAIGLDKDKP  NRVTKVAVKM  LKSDATEKDL
301  SDLISEMEM  M  KMIGKHKNII  NLLGACTQDG  PLYVIMEYAS  KGNLREYLQA  RPPGLEICY
361  NP SHNPEEQ  L  SSKDLVSCAY  QVARGMEYLA  SKKCIHRDLA  ARNVLVTE  DN  VMKIADFGLA
421  RDIHHIDY  YK  KTTNGRLPVK  WMAPEALFDR  IYTHQSDVWS  FGVLLWEIF  T  LGGSPYPGPV
481  VEELFKLL  KE  GHRMDKPSNC  TNELYMMMRD  CWHAVPSQRP  TFKQLVEDLD  RIVALTSNQE
  
```

#### Recombinant FGFR1 (V561M) nucleotide sequence:

```

1  atgtccccta  tactaggtta  ttggaaaatt  aagggccttg  tgcaaccac  tcgacttctt
61  ttggaatata  ttgaagaaaa  atatgaagag  catttgtag  agcgcgatga  aggtgataaa
121  tggcgaagaa  aaaagtttga  attgggtttg  gagtttcca  atcttcctta  ttatatgtat
181  ggtgatgtta  aattaacaca  gtctatggcc  atcatacgtt  atatagctga  caagcacaac
241  atgttgggtg  gttgtccaaa  agagcgtgca  gagatttcaa  tgcttgaag  agcggttttg
301  gatattagat  acggtgtttc  gagaattgca  tatagtaaag  actttgaaac  tctcaaagtt
361  gattttctta  gcaagctacc  tgaaatgctg  aaaatgttcg  aagatcgttt  atgtcataaa
421  acataattaa  atggtgatca  tgtaacccat  cctgacttca  tgttgatga  cgctcttgat
481  gtgttttat  acatggacc  aatgtgcctg  gatgcgttcc  caaaattagt  ttgttttaa
541  aaacgtattg  aagctatccc  acaaattgat  aagtacttga  aatccagcaa  gtatatagca
601  tggcctttgc  agggctggca  agccacgttt  ggtggtggcg  accatcctcc  aaaatcggat
661  ctggaagtcc  tgttccaggg  gcccgaaatt  atgctagcag  gggctctctga  gtatgagctt
721  cccgaagacc  ctgcctggga  gctgcctcgg  gacagactgg  tcttaggcaa  accctggga
781  gagggctgct  ttgggcagg  ggtggtggca  gaggctatcg  ggctggaca  agcacaaccc
841  aaccgtgtga  ccaaagtggc  tgtgaagatg  ttgaagtcgg  acgcaacaga  gaaagacttg
901  tcagacctga  tctcagaaat  ggagatgatg  aagatgatcg  ggaagcataa  gaatatcatc
961  aacctgctgg  gggcctgcac  gcaggatggt  cccttgtag  tcatcatgga  gtatgcctcc
1021  aagggcaacc  tgcgggagta  cctgcaggcc  cggaggccc  cagggctgga  atactgctac
1081  aacccagacc  acaaccaga  ggagcagctc  tctccaagg  acctggtgtc  ctgcccctac
1141  caggtggccc  gaggcattga  gtatctggcc  tccaagaagt  gcatacaccg  agacctggca
1201  gccaggaatg  tcctggtgac  agaggacaat  gtgatgaaga  tagcagactt  tggcctcgca
1261  cgggacattc  accacatcga  ctactataaa  aagacaacca  acggccgact  gctgtgaag
1321  tggatggcac  cggaggcatt  atttgaccgg  atctacacc  accagagtga  tgtgtgtct
1381  ttcgggggtg  tcctgtggga  gatcttcaact  ctggggcggc  ccccatacc  cgggtgtgct
1441  gtggaggaac  ttttcaagct  gctgaaggag  ggtcacccga  tggacaagcc  cagtaactgc
1501  accaacgagc  tgtacatgat  gatgcgggac  tgctggcatg  cagtgcctc  acagagacc
1561  acctcaagc  agctggtgga  agacctggac  cgcatcgtgg  ccttgacct  caaccaggag
1621  taa
  
```

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.