

## Certificate of Analysis

### EGFR (L858R), active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-626, 14-626-K, 14-626M

Parent Lot # 31001U

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 EGFR, amino acids 696–end containing the mutation L858R. This mutation has been associated with non-small cell lung cancer patients who demonstrate clinical responsiveness to the tyrosine kinase inhibitor gefitinib (Iressa, ZD1839). Expressed by baculovirus in Sf21 insect cells. Purified using glutathione-agarose. Purity 73% by SDS-PAGE and Coomassie blue staining. MW = 85.8kDa.

**Specific Activity (Parent lot# 31001U):** 376U/mg, where one unit of EGFR activity is defined as 1nmol phosphate incorporated into 0.1mg/ml poly(Glu, Tyr) (4:1) per minute at 30°C with a final ATP concentration of 100µM.

**Formulation:** 2.2mg/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.

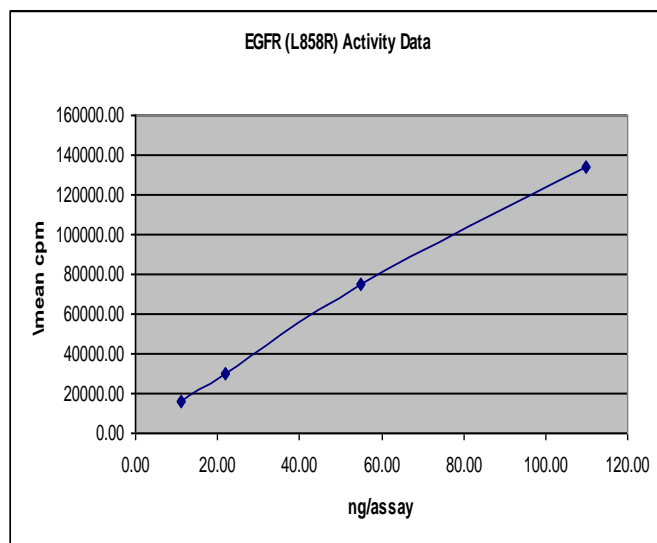
**Storage and Stability:** On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 6 months from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the 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.

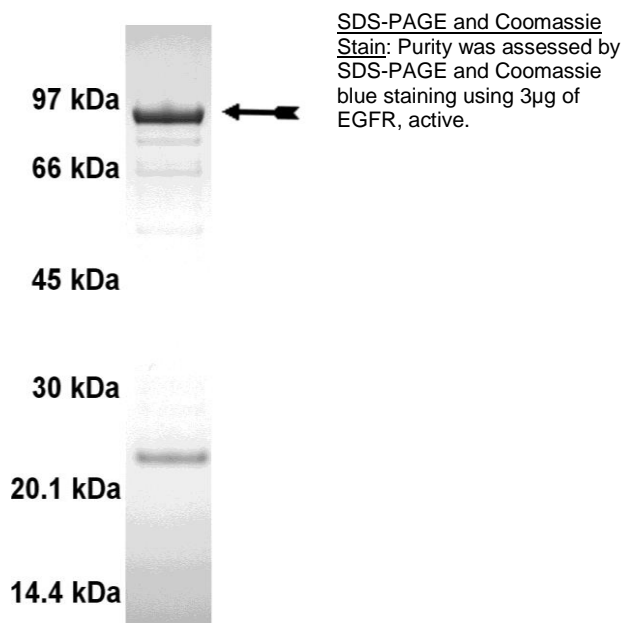
**FOR IN VITRO RESEARCH USE ONLY  
NOT FOR USE IN HUMANS OR ANIMALS**

### Quality Control Testing

**Kinase Assay:** 11–110ng of this lot of enzyme phosphorylated 0.1mg/ml poly(Glu,Tyr) (4:1) 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 identity as EGFR 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. **Poly(Glu, Tyr) (4:1):** Use at a final assay concentration of 0.1mg/ml. Prepare a 1mg/ml stock. Add 2.5µl of stock per assay point.
3. **EGFR, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 11–110ng 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 5µl of 10 x reaction buffer per assay to wells
2. Add 2.5µl of **poly(Glu, Tyr) (4:1)**.
3. Add **2.5µl (11–110ng) EGFR, active**.
4. Add 5µl of dH<sub>2</sub>O.
5. Add 10µl of diluted [ $\gamma$ -<sup>33</sup>P]ATP mixture.
6. Incubate for 10 minutes at 30°C.
7. Stop the reaction by adding 5µl of 3% phosphoric acid.
8. Transfer a 10µl aliquot onto the appropriate area of a **Filtermat A**.
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 1µl of 30% phosphoric acid.

## Certificate of Analysis

### EGFR (L858R) Sequence Information

<b><u>Protein</u></b>	Human EGFR (L858R)
<b><u>Tags</u></b>	N-Terminal GST
<b><u>Native sequence</u></b>	G241 of recombinant sequence is equivalent to G696 of native human EGFR
<b><u>Accession number</u></b>	GenBank X00588

#### Recombinant EGFR amino acid sequence:

```

1  MSPILGYWKI  KGLVQPTRL  LEYLEEKYEE  HLYERDEGDK  WRNKKFELGL  EFPNLPYYID
61  GDVKLTQ SMA  IIRYIADKHN  MLGGCPKERA  EISMLEGAVL  DIRYGVSRIA  YSKDFETLKV
121  DFLSKLP EML  KMFEDRLCHK  TYLNGDHVTH  PDFMLYDALD  VVLYMDPMCL  DAFPKLVCFK
181  KRIEAI PQID  KYLKSSKYIA  WPLQGWQATF  GGGDHPPKSD  LEVLFQGP EF  KGMGIRNSKG
241  GEAPNQALLR  ILKETEFK KI  KVLGSGAFGT  VYKGLWIPEG  EKVKIPVAIK  ELREATSPKA
301  NKEILDEAYV  MASVDNPHVC  RLLGICLTST  VQLITQLMPF  GCLLDYVREH  KDNIGSQYLL
361  NWCVQIAKGM  NYLED RRLVH  RDLAARNVLV  KTPQHVKITD  FGRAKLLGAE  EKEYHAEGGK
421  VPIKWMALES  ILHRIYTHQS  DVWSYGVTVW  ELMTFGSKPY  DGIPASEISS  ILEKGERLPQ
481  PPICTIDVYM  IMVKCWMIDA  DSRPKFRELI  IEF SKMARDP  QRYLVIQ GDE  RMHLPSP TDS
541  NFYRALMDEE  DMDDV DDADE  YLIPQQGFFS  SPSTSRTPLL  SLSATSNN S  TVACIDR NGL
601  QSCPIKEDSF  LQRYSSDPTG  ALTEDSIDDT  FLPVPEYINQ  SVPKRPAGSV  QNPVYHNQPL
661  NPAPSRDPHY  QDPHSTAVGN  PEYLN TVQPT  CVNSTFDSPA  HWAQKGS HQI  SLDNPDYQQD
721  FFPKEAKPNG  IFKGSTAENA  EYLRVAPQSS  EFIGA

```

#### Recombinant EGFR nucleotide sequence:

```

1  atgtccccta  tactagg tta  ttggaaaatt  aagggccttg  tgcaaccac  tcgacttctt
61  ttggaatata  ttgaagaaa  atatgaagag  catttgtatg  agcgcgatga  aggtgataaa
121  tggcgaaaca  aaaagtttga  attgggtttg  gagtttccca  atcttcctta  ttatattgat
181  ggtgatg tta  aattaacaca  gtctatggcc  atcatacggt  atatagctga  caagcacaac
241  atgttgggtg  gttgtccaaa  agagcgtgca  gagatttcaa  tgcttgaagg  agcggttttg
301  gatattagat  acggtgtttc  gagaattgca  tatagtaaag  actttgaaac  tctcaaagtt
361  gattttctta  gcaagctacc  tgaaatgctg  aaaatgttcg  aagatcgttt  atgtcataaa
421  acatatttaa  atggtgatca  tgtaaccat  cctgacttca  tgttgtatga  cgctcttgat
481  gttgttttat  acatggacc  aatgtgectg  gatgcgttcc  caaaattagt  ttgttttaa
541  aaacgtattg  aagctatccc  acaaattgat  aagtacttga  aatccagcaa  gtatatagca
601  tggcctttgc  agggctggca  agccacgttt  ggtggtggcg  accatcctcc  aaaatcggat
661  ctggaagtgc  tgttccagg  gcccgattc  aaaggcatgg  ggatccggaa  ttcaaagggt
721  ggagaagctc  ccaaccaagc  tctcttgagg  atcttgaagg  aaactgaatt  caaaaagatc
781  aaagtgtctg  gctccggtgc  gttcggcagc  gtgtataagg  gactctggat  cccagaaggt
841  gagaaagtta  aaattcccgt  cgctatcaag  gaattaagag  aagcaacatc  tccgaaagcc
901  aacaaggaaa  tcctcgatga  agcctacgtg  atggccagcg  tggacaacc  ccagtggtgc
961  cgctgctg  gcatctgcct  cacctccacc  gtgcaactca  tcacgcagct  catgcccttc
1021  ggetgcctcc  tggactatgt  ccgggaacac  aaagacaata  ttggctccca  gtacctgctc
1081  aactggtgtg  tgcagatcgc  aaagggcatg  aactacttgg  aggaccgtcg  cttggtgcac
1141  cgcgacctgg  cagccaggaa  cgtactggtg  aaaacaccgc  agcatgtcaa  gatcacagat
1201  tttgggctgg  ccaaactgct  ggggtcggaa  gagaaagaat  accatgcaga  aggaggcaaa
1261  gtgcctatca  agtggtatgg  attggaatca  attttacaca  gaatctatac  ccaccagagt
1321  gatgtctgga  gctacggggt  gaccgtttgg  gagttgatga  ctttggatc  caagccatat
1381  gacggaatcc  ctgccagcga  gatctcctcc  atcctggaga  aaggagaacg  cctccctcag
1441  ccaccatata  gtaccatcga  tgtctacatg  atcatggtca  agtgctggat  gatagacgca
1501  gatagtgc  caaagttccg  tgagttgatc  atcgaattct  caaaatggc  ccgagacccc
1561  cagcgtacc  ttgtcattca  gggggatgaa  agaatgcatt  tgccaagtcc  tacagactcc
1621  aacttctacc  gtgcctgat  ggatgaagaa  gacatggacg  acgtggtgga  tgccgacgag
1681  tacctcatcc  cacagcagg  cttcttcagc  agcccctcca  cgtcacggac  tcccctcctg

```

## Certificate of Analysis

```
1741 agctctctga gtgcaaccag caacaattcc accgtggctt gcattgatag aaatgggctg
1801 caaagctgtc ccatcaagga agacagcttc ttgcagcgat acagctcaga cccacagggc
1861 gccttgactg aggacagcat agacgacacc ttctctccag tgcctgaata cataaaccag
1921 tccgttccca aaaggcccgc tggtctctgt cagaatcctg tctatcacia tcagcctctg
1981 aaccccgcgc ccagcagaga cccacactac caggaccccc acagcactgc agtgggcaac
2041 cccgagtatc tcaacactgt ccagcccacc tgtgtcaaca gcacattcga cagccctgcc
2101 cactgggccc agaaaggcag ccaccaaatt agcctggaca accctgacta ccagcaggac
2161 ttctttccca aggaagccaa gccaaatggc atctttaagg gctccacagc tgaaaatgca
2221 gaatacctaa ggtcgcgcc acaaagcagt gaatttattg gagcatga
```

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.