

Certificate of Analysis

EGFR (T790M,L858R), active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-721, 14-721-K, 14-721M

Parent Lot # 1691467

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 mutations T790M and L858R, expressed by baculovirus in Sf21 insect cells. Purified using glutathione-agarose. Purity 81.9% by SDS-PAGE and Coomassie blue staining. MW = 85.8kDa.

Specific Activity (Parent lot# 1691467): 107U/mg, where one unit of EGFR (T790M,L858R), active activity is defined as 1nmol phosphate incorporated into 250 μ M (GGMEDIYFEFMGGKKK) per minute at 30°C with a final ATP concentration of 100 μ M.

Formulation: 0.501mg/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.

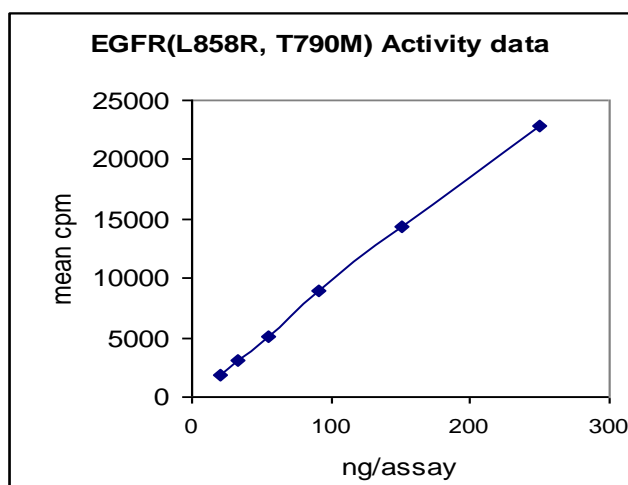
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

Handling Recommendations: Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled 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: 20–251ng of this lot of enzyme phosphorylated 250 μ M (GGMEDIYFEFMGGKKK) 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 (T790M, L858R), 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. (GGMEDIYFEFMGGKKK):** Use at a final assay concentration of 250 μ M. Make up a 2.5mM stock. Add 2.5 μ l of stock per assay point.
- 3. EGFR (T790M, L858R), active:** Dilute with 20mM MOPS-NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 20–251 ng 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:

1. Add 5 μ l of 5 x reaction buffer per assay to wells.
2. Add 2.5 μ l of (GGMEDIYFEFMGGKKK).
3. Add **2.5 μ l (20–251ng) EGFR (T790M, L858R) 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 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 μ l of 30% phosphoric acid.

Certificate of Analysis

EGFR-(696–end, T790M, L858R) Sequence Information

Protein Human EGFR (T790M,L858R)

Tags N-terminal GST

Native sequence G241 of the recombinant protein is equivalent to G696 of human EGFR (T790M,L858R). The amino acid substitution L858R is one of several heterozygous mutations that have been identified in Non-Small-Cell Lung Cancer (NSCLC) patients who have clinical responses to the EGFR inhibitor Iressa®. There is some evidence that these mutations result in elevated activity and enhanced sensitivity to Iressa® (Lynch *et al.*, (2004), N. Engl. J. Med. **350**, 2129-2139). In patients with tumours bearing Iressa®-sensitive mutations, resistant subclones containing an additional EGFR mutation, T790M, emerge in the presence of the drug (Pao *et al.*, (2005), PLoS Medicine **2**, 225-235; Kwak *et al.*, (2005), Proc. Natl. Acad. Sci. USA **102**, 7665-7670.). It has been shown experimentally that the T790M mutation leads to high-level functional resistance to Iressa® (Kobayashi *et al.*, (2005), Cancer Res. **65**, 7096-7101.)

Accession number GenBank X00588

Recombinant EGFR (T790M, L858R) amino acid sequence:

```

1 MSPILGYWKI KGLVQPTRLL LEYLEEKYEE HLYERDEGDK WRNKKFELGL EFPNLPYYID
61 GDVKLTQSM A IIRYIADKHN MLGGCPKERA EISMLEGAVL DIRYGVSR IA YSKDFETLKV
121 DFLSKLPEML KMFEDRLCHK TYLNGDHVTH PDFMLYDALD VVLYMDPMCL DAFPKLVCFK
181 KRIEAI PQID KYLKSSKYIA WPLQGWQATF GGGDHPPKSD LEVLFQGP EF KGMGIRNSKG
241 GEAPNQALLR ILKETEFKKI KVLGSGAFGT VYKGLWIPEG EKVKIPVAIK ELREATSPKA
301 NKEILDEAYV MASVDNPHVC RLLGICLTST VQLIMQLMPF GCLLDYVREH KDNIGSQYLL
361 NWCVQIAKGM NYLED RRLVH RDLAARNVLV KTPQHVKITD FGRAKLLGAE EKEYHAEGGK
421 VPIKWMAL ES ILHRIYTHQS DVWSYGVTVW ELMTFGSKPY DGIPASEISS ILEKGERLPQ
481 PPICTIDVYM IMVKCWMIDA DSRPKFRELI IEFKSMARDP QRYLVIQGDE RMHLPSP TDS
541 NFYRALMDEE DMDDVVD ADE YLIPQQGFFS SPSTSRTPLL SLSATSNN S TVACIDRNG L
601 QSCPIKEDSF LQRYSSDPTG ALTEDSIDDT FLPVPEYINQ SVPKRPAGSV QNPVYHNQPL
661 NPAPSRDPHY QDPHSTAVGN PEYLN TVQPT CVNSTFDSPA HWAQK GSHQI SLDNPDYQQD
721 FFPKEAKPNG IFKGSTAENA EYLRVAPQSS EFIGA
  
```

Recombinant EGFR (T790M, L858R) nucleotide sequence:

```

1 atgtccccta tactaggtta ttgaaaatt aagggccttg tgcaaccac tcgacttctt
61 ttggaatata ttgaagaaaa atatgaagag catttगतatg agcgcgatga aggtgataaa
121 tggcgaacaa aaaagtttga attgggtttg gagtttcca atcttcctta ttatattgat
181 ggtgatgtta aattaacaca gtctatggcc atcatacgtt atatagctga 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 tgttगतatga cgctcttgat
481 gttgttttat acatggacc aatgtgcctg gatgcgttcc caaaattagt ttgttttaa
541 aaactgattg aagctatccc acaaattgat aagtacttga aatccagcaa gtatatagca
601 tggcctttgc agggctggca agccacgtt ggtgggtggcg accatcctcc aaaatcggat
661 ctggaagttc tgttccagg gcccgaattc aaaggcatgg ggatccggaa ttcaaagggt
721 ggagaagctc ccaaccaagc tctcttgagg atcttgaagg aaactgaatt caaaaagatc
781 aaagtगतtg gctccggtgc gttcggcacg gtgtataagg gactctggat cccagaaggt
841 gagaaagtta aaattcccgt cgctatcaag gaattaagag aagcaacatc tccgaaagcc
901 aacaaggaaa tcctcगतga agcctacgtg atggccagcg tggacaacc ccacgtगतc
961 gcctगतtg gcatctgcct cacctcacc gtgcaactca tcatgcagct catgccctt
  
```

Certificate of Analysis

```
1021 ggctgcctcc tggactatgt ccgggaacac aaagacaata ttggctccca gtacctgctc
1081 aactggtgtg tgcagatcgc aaagggcatg aactacttgg aggaccgtcg cttggtgcac
1141 cgcgacctgg cagccaggaa cgtactggtg aaaacaccgc agcatgtcaa gatcacagat
1201 tttgggcggg ccaaactgct ggggtcggaa gagaaagaat accatgcaga aggaggcaaa
1261 gtgcctatca agtggatggc atttgaatca attttacaca gaatctatac ccaccagagt
1321 gatgtctgga gctacggggg gaccgttgg gagttgatga cctttggatc caagccatat
1381 gacggaatcc ctgccagcga gatctcctcc atcctggaga aaggagaacg cctccctcag
1441 ccacccatat gtaccatcga tgtctacatg atcatggtca agtgctggat gatagacgca
1501 gatagtcgcc caaagttccg tgagttgatc atcgaattct ccaaaatggc ccgagacccc
1561 cagcgtacc ttgtcattca gggggatgaa agaatgcatt tgccaagtcc tacagactcc
1621 aacttctacc gtgccctgat ggatgaagaa gacatggacg acgtggtgga tgccgacgag
1681 tacctcatcc cacagcaggg cttcttcagc agcccctcca cgtcacggac tcccctcctg
1741 agctctctga gtgcaaccag caacaattcc accgtggctt gcattgatag aaatgggctg
1801 caaagctgtc ccatcaagga agacagctt ttgcagcga acagctcaga cccacaggc
1861 gccttgactg aggacagcat agacgacacc ttctcccag tgcctgaata cataaaccag
1921 tccgttccca aaaggccgc tggctctgtg cagaatcctg tctatcacia tcagcctctg
1981 aaccccgccg 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 gggtcgcgcc 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