

## Certificate of Analysis

### PDGFR alpha (550-end), active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-467, 14-467-K, 14-467M

Parent Lot # 1767085

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 6His-tagged recombinant human PDGFR alpha residues 550–end. Expressed by baculovirus in Sf21 insect cells. Purified using Ni<sup>2+</sup>/NTA-agarose. Purity 80.3% by SDS-PAGE and Coomassie blue staining. MW = 63.5kDa.

**Specific Activity (Parent lot# 1767085):** 29U/mg, where one unit of PDGFR alpha, active 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:** 1.127mg/ml of enzyme in 50mM Tris/HCl pH7.5, 0.1mM EGTA, 150mM NaCl, 0.03% Brij-35, 270mM sucrose, 0.2mM PMSF, 1mM benzamidine, 0.1% 2-mercaptoethanol. Frozen solution.

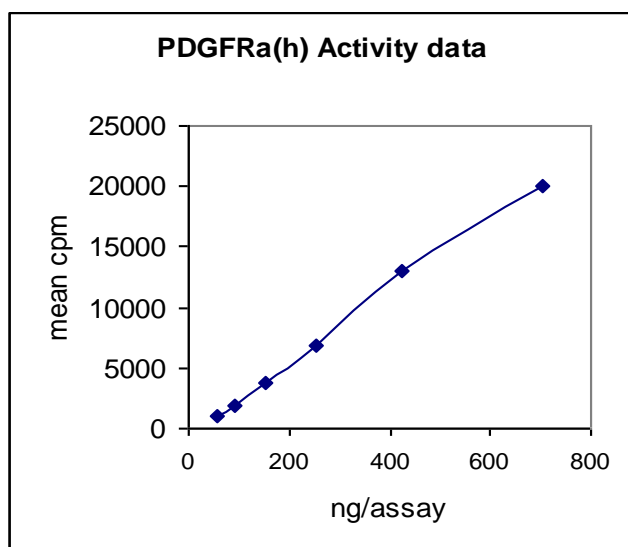
**Storage and Stability:** On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

**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**

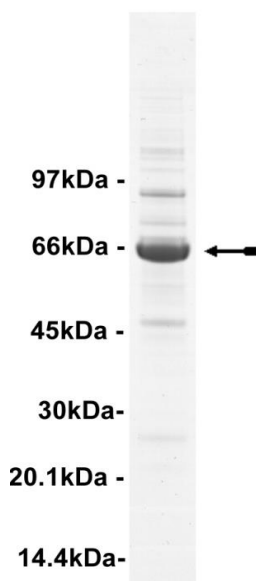
### Quality Control Testing

**Kinase Assay:** 55–704ng of this lot of enzyme phosphorylated 0.1mg/ml poly(Glu, Tyr) 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 PDGFR alpha with the translated sequence listed on page three.

**SDS-PAGE and Coomassie Stain:** Purity was assessed by SDS-PAGE and Coomassie blue staining using 3µg of PDGFR alpha, active.



## Certificate of Analysis

### Kinase Assay Protocol

#### Stock Solutions:

1. **5 x Reaction Buffer:** 40mM MOPS/NaOH pH7.0, 1mM EDTA.
2. **MnCl<sub>2</sub>:** Use at a final assay concentration of 10mM. Prepare a 200mM stock in dH<sub>2</sub>O and add 1.25µl of stock per assay point.
3. **Poly(Glu, Tyr) (4:1):** Use at a final assay concentration of 0.1mg/ml. Prepare a 1mg/ml stock and add 2.5µl of stock per assay point.
4. **PDGFR alpha, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1 % 2-mercaptoethanol, 1mg/ml BSA. Use 55–704ng per assay point.
5. **[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 5µl of 5 x reaction buffer per assay to wells.
2. Add 2.5µl of **poly(Glu, Tyr) (4:1)**.
3. Add 1.25µl of MnCl<sub>2</sub>.
4. Add **2.5µl (55–704ng) PDGFR alpha, active**.
5. Add 3.75µl of dH<sub>2</sub>O.
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 of 3% phosphoric acid.
9. Transfer a 10µl aliquot onto the appropriate area of a **Filtermat A**.
10. Wash the filtermat three times for 5 minutes with 75mM phosphoric acid.
11. Wash the filtermat once for 2 minutes with methanol.
12. 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

## PDGFR alpha Sequence Information

<b>Protein</b>	human PDGFR alpha
<b>Tags</b>	N-terminal 6His
<b>Native sequence</b>	K16 of the recombinant protein is equivalent to K550 of human PDGFR alpha
<b>Accession number</b>	GenBank M21574

### Recombinant PDGFR alpha amino acid sequence:

```

1 MAHHHHHHEN LYFQ GKQKPR YEIRWRVIES ISPDGHEIY VDPMLPYDS RWEFPRDGLV
61 LGRVLGSGAF GKVEGTAYG LRSQPVMKV AVKMLKPTAR SSEKQALMSE LKIMTHLGPH
121 LNIIVLLGAC TKS GPIYIIT EYCFYGLLVN YLHKNRDSFL SHHPEKPKKE LDIFGLNPAD
181 ESTRSYVILS FENNGDYMDM KQADTTQYVP MLERKEVSKY SDIQRSLYDR PASYKKKSML
241 DSEVKNLLSD DNSEGLTLLD LLSFTYQVAR GMEFLASKNC VHRDLAARNV LLAQGKIVKI
301 CDFGLARDIM HDSNYVSKGS TFLPVKMAP ESIFDNLYTT LSDVWSYGIL LWEIFSLGGT
361 PYPGMMVDST FYNKIKSGYR MAKPDHATSE VYEIMVKCWN SEPEKRPSFY HLSEIVENLL
421 PGQYKKSIEK IHLDFLKSDH PAVARMRVDS DNAYIGVTYK NEEDKLDWE GGLDEQRSLA
481 DSGYIIPDPD IDPVEEEDL GKRNRHSSQT SEESAIETGS SSSTFIKRED ETIEDIDMMD
541 DIGIDSSDLV EDSFL
  
```

### Recombinant PDGFR alpha nucleotide sequence:

```

1 atggcgcatt accatcacca tcatgaaaac ctgtattttc agggcaaaca gaaaccgagg
61 tatgaaattc gctggagggt cattgaatca atcagcccgg atggacatga atatatttat
121 gtggaccgga tgcagctgcc ttatgactca agatgggagt ttccaagaga tggactagtg
181 cttggctggg tcttggggctc tggagcgttt gggaaaggagg ttgaaggaac agcctatgga
241 ttaagccggt cccaacctgt catgaaagt gcaagtgaaga tgctaaaacc cacggccaga
301 tccagtgaaa aacaagctct catgtctgaa ctgaagataa tgactcacct ggggccacat
361 ttgaacattg taaacttctt gggagcctgc accaagtcag gccccattta catcatcaca
421 gagtattgct tctatggaga tttggtaaac tttttgcata agaataggga tagcttcctg
481 agccaccacc cagagaagcc aaagaaagag ctggatatct ttggattgaa ccctgctgat
541 gaaagcacac ggagctatgt tattttatct tttgaaaaca atggtgacta catggacatg
601 aagcaggctg atactacaca gtatgtccc atgctagaaa ggaaagagggt ttctaaatat
661 tccgacatcc agagatcact ctatgatcgt ccagcctcat ataagaagaa atctatgtta
721 gactcagaag tcaaaaacct ctttcagat gataactcag aaggccttac tttattggat
781 ttgttgagct tcacatatca agttgccga ggaatggagt ttttggcttc aaaaaattgt
841 gtcaccgctg atctggctgc tcgcaacgct ctcctggcac aaggaaaaat tgtgaagatc
901 tgtgactttg gcctggccag agacatcatg catgattcga actatgtgtc gaaaggcagt
961 acctttctgc ccgtgaagtg gatggctcct gagagcatct ttgacaacct ctacaccaca
1021 ctgagtgatg tctgttctta tggcattctg ctctgggaga tcttttccct tgggtggacc
1081 ccttaccgag gcatgatggg ggattctact ttctacaata agatcaagag tgggtaccgg
1141 atggccaagc ctgaccacgc taccagtga gctctacgaga tcatgggtgaa atgctggaac
1201 agtgagccgg agaagagacc ctccttttac cacctgagtg agattgtgga gaatctgctg
1261 cctggacaat ataaaaagag ttatgaaaaa attcacctgg acttctgtaa gaggaccat
1321 cctgctgtgg cacgcatgct tgtggactca gacaatgcat acattgggtg cacctacaaa
1381 aacgaggaag acaagctgaa ggactgggag ggtggctctg atgagcagag actgagcgtc
1441 gacagtggct acatcattcc tctgcctgac attgaccctg tccctgagga ggaggacctg
1501 ggcaagagga acagacacag ctgcagacc tctgaagaga gtgccattga gacgggttcc
1561 agcagttcca cttcatcaa gagagaggac gagaccattg aagacatcga catgatggac
1621 gacatcggca tagactcttc agacctggtg gaagacagct tcctgtaa
  
```

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.