

Certificate of Analysis

HePTP, active

(Recombinant enzyme expressed in *E. coli* cells)

Item # 14-593, 14-593-K, 14-593M

Parent Lot # 1922202

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 HePTP, amino acids 22–end, expressed in *E. coli* cells. Purified using Ni²⁺/NTA agarose. Purity 72.9% by SDS-PAGE and Coomassie blue staining. MW = 40.7kDa.

Specific Activity (Parent lot# 1922202): 8316U/mg, where one unit of HePTP, active activity is defined as the release of 1nmol of phosphate per minute from the phosphorylated substrate 6,8-difluoro-4-methylumbelliferyl phosphate (DiFMUP) at room temperature.

Formulation: 0.812mg/ml of enzyme in 50mM Tris/HCl pH7.5, 150mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 10% glycerol, 0.2mM PMSF, 1mM benzamidine, 0.1% 2-mercaptoethanol. Frozen solution.

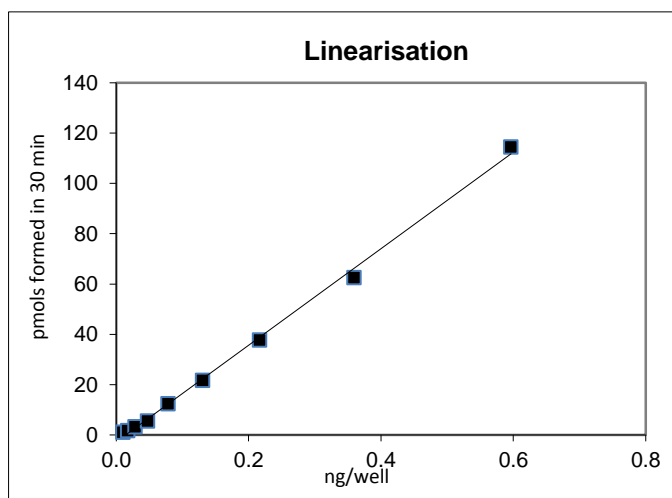
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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.

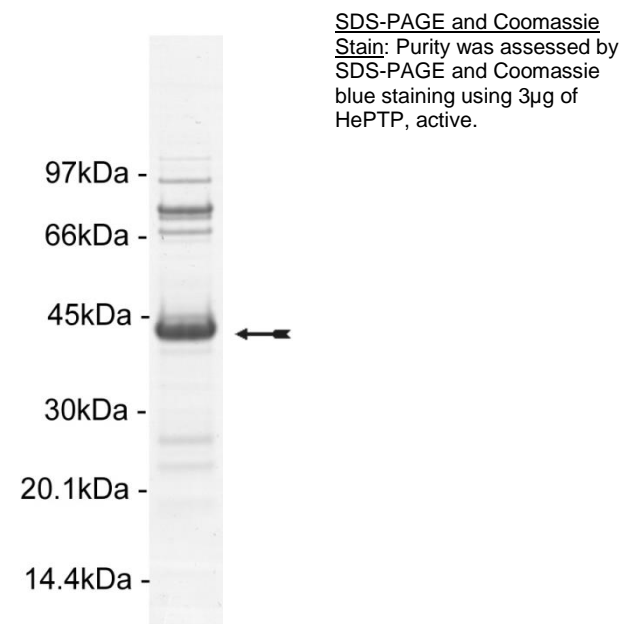
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Phosphatase Assay: 0.01–0.6ng of this lot of enzyme dephosphorylated 200µM DiFMUP in the assay described on page two. Assay background was subtracted from the actual Fluorescence Intensity (FI) to yield the results shown below. Quantification of FI was against a 6,8-difluoro-7-hydroxy-4-methylcoumarin (DiFMU) standard curve.



MS Tryptic Fingerprint: Confirmed identity as HePTP with the translated sequence listed on page three.



Certificate of Analysis

Phosphatase Assay Protocol

Stock Solutions:

1. **Reaction Buffer:** 60mM Hepes pH7.2, 150mM NaCl, 1mM EDTA, 0.17mM DTT, 0.83 (v/v)% glycerol, 0.017 (w/v)% BSA, 0.002% Brij-35.
2. 500 μ M DiFMUP (Molecular Probes Catalogue# D6567) in water.
3. 100mM sodium orthovanadate.
4. 500 μ M DiFMU (Molecular Probes Catalogue# D6566) in water for the calibration curve.

Assay Procedure:

1. Dilute **HePTP** in reaction buffer and use 0.01–0.6ng in 15 μ l per assay point.
2. Add 10 μ l DiFMUP 500 μ M stock solution (200 μ M final assay concentration).
3. Incubate for 30 minutes at room temperature.
4. Stop the reaction by adding 5 μ l of 100mM sodium orthovanadate.
5. Read FI using an appropriate reader (Excitation 340nm; Emission 450nm).
6. Subtract the zero enzyme values from each FI reading and calculate the enzyme activity by conversion to nmoles product formed using a DiFMU standard calibration curve.

Certificate of Analysis

HePTP Sequence Information

<u>Protein</u>	HePTP
<u>Tags</u>	N-terminal 6His
<u>Native sequence</u>	M21 of the recombinant protein is equivalent to M22 of human HePTP
<u>Accession number</u>	GenBank NM_002832

Recombinant HePTP amino acid sequence:

```

1  MGSSHHHHHH  SSGLVPRGSH  MTQPPPEKTP  AKKHVRLQER  RGSNVALMLD  VRSLGAVEPI
61  CSVNTPREVT  LHFLRTAGHP  LTRWALQRQP  PSPKQLEEEF  LKIPSNFVSP  EDLDIPGHAS
121  KDRYKTIILPN  PQSRVCLGRA  QSQEDGDYIN  ANYIRGYDVK  EKVIYIATQGP  MPNTVSDSDFWE
181  MVWQEEVSLI  VMLTQLREGK  EKCIVHYWPT  EETYGPFQIR  IQDMKECPEY  TVRQLTIQYQ
241  EERRSVKHIL  FSAWPDHQT  ESAGPLLRLV  AEVEESPETA  AHPGPIVVHC  SAGIGRTGCF
301  IATRIGCQQL  KARGEVDILG  IVCQLRLDRG  GMIQTAEQYQ  FLHHTLALYA  GQLPEEPS
  
```

Recombinant HePTP nucleotide sequence:

```

1  atgggcagca  gccatcatca  tcatcatcac  agcagcggcc  tgggtgccg  cggcagccat
61  atgaccagc  ctccgcctga  aaaaacgcca  gccagaagc  atgtgcgact  gcaggagag
121  cggggctcca  atgtgctct  gatgctggac  gttcggctcc  tgggggccc  agaaccatc
181  tgctctgtga  acacacccc  ggaggtcacc  ctacactttc  tgcgcactgc  tggacacccc
241  ctaccgcgt  gggcccttca  gcgccagcca  cccagcccca  agcaactgga  agaagaattc
301  ttgaagatcc  cttcaaactt  tgtcagcccc  gaagacctgg  acatccctgg  ccacgcctcc
361  aaggaccgat  acaagaccat  cttgccaaat  ccccagagcc  gtgtctgtct  aggccgggca
421  cagagccagg  aggacggaga  ttacatcaat  gccaaactaca  tccgaggcta  tgacgggaag
481  gagaaggtct  acattgccac  ccagggcccc  atgcccaaca  ctgtgtcgga  cttctgggag
541  atggtgtggc  aagaggaagt  gtccctcatt  gtcatgctca  ctcagctccg  agagggcaag
601  gagaaatgtg  tccactactg  gccacagaa  gaggaacct  atggaccctt  ccagatccgc
661  atccaggaca  tgaagagtg  cccagaatac  actgtgcggc  agctcaccat  ccagtaccag
721  gaagagcgcc  ggtcagtaaa  gcacatcctc  ttttcggcct  ggccagacca  tcagacacca
781  gaatcagctg  ggcccctgct  gcgcctagt  gcagaggtgg  aggagagccc  ggagacagcc
841  gccacccc  ggccctatcgt  agtccactgc  agtgcaggg  ttggccggac  gggctgcttc
901  atcgccagc  gaattggctg  tcaacagctg  aaagcccag  gagaagtgga  cattctgggt
961  attgtgtgcc  aactgcccgt  agacagaggg  gggatgatcc  agacggcaga  gcagtaccag
1021  ttctgcacc  acactttggc  cctgtatgca  ggccagctgc  ctgaggaacc  cagcccctga
  
```

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.