

Certificate of Analysis

DUSP22, active

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

Item # 14-641, 14-641-K, 14-641M

Parent Lot # 1922204

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 DUSP22, full length, expressed in *E. coli* cells. Purified using glutathione agarose. Purity 91.9% by SDS-PAGE and Coomassie blue staining. MW = 47.3kDa.

Specific Activity (Parent lot# 1922204): 360U/mg, where one unit of DUSP22, 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: 1.629mg/ml of enzyme in 50mM Tris/HCl pH7.5, 150mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 1mM benzamidine, 0.2mM PMSF, 0.1% 2-mercaptoethanol. Frozen solution.

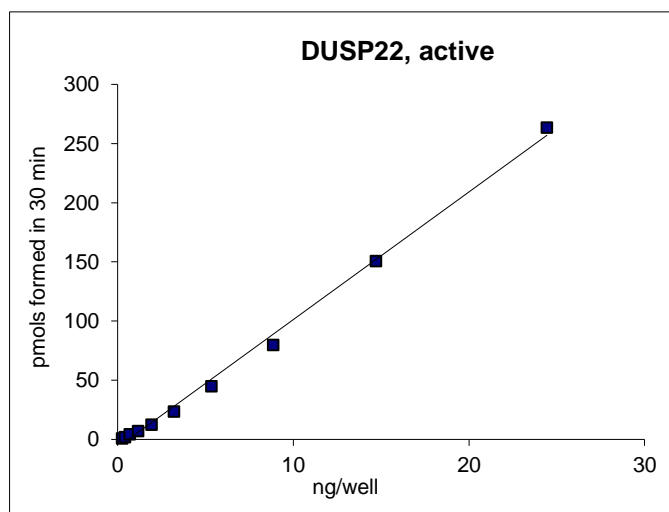
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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.

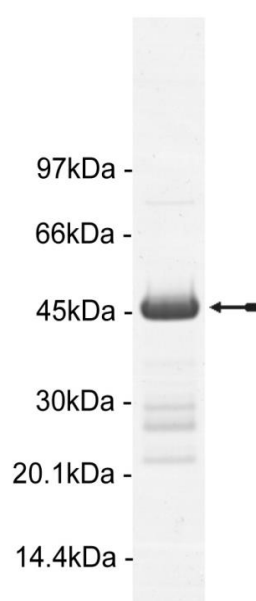
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Phosphatase Assay: 0.2–2.4ng 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 DUSP22 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 DUSP22, active.

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. 500µM DiFMUP (Molecular Probes Catalogue# D6567) in water.
3. 100mM sodium orthovanadate.
4. 500µM DiFMU (Molecular Probes Catalogue# D6566) in water for the calibration curve.

Assay Procedure:

1. Dilute **DUSP22** in reaction buffer and use 0.2–2.4ng in 15µl per assay point.
2. Add 10µl DiFMUP 500 µM stock solution (200µM final assay concentration).
3. Incubate for 30 minutes at room temperature.
4. Stop the reaction by adding 5µl 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.

Preparation of DiFMU Standard Calibration Curve

1. Prepare a series of standards according to the table below:

Volume of 100µM DiFMU (product)	Volume of 100µM DiFMUP (substrate)	pmoles DiFMU per well	Represents Substrate Conversion of:
0µl	100µl	0	0%
1µl	99µl	25	1%
2µl	98µl	50	2%
3µl	97µl	75	3%
4µl	96µl	100	4%
5µl	95µl	125	5%
8µl	92µl	200	8%
12µl	88µl	300	12%

2. Add 25µl of each prepared standard to triplicate wells and read fluorescence intensity using an appropriate reader (Excitation 358nm, Emission 455nm).

Note: Substrate conversion above 15% introduces the possibility of skewing the curve so that calculated specific activity will be less than actual specific activity. This substrate limitation kinetic varies with each enzyme and may be determined by plotting product formed per unit time versus enzyme concentration. At low enzyme concentrations, the rate of product formation will be proportional to the enzyme concentration used. This relationship plateaus at higher enzyme concentrations.

Certificate of Analysis

DUSP22 Sequence Information

<u>Protein</u>	Human DUSP22
<u>Tags</u>	N-Terminal GST
<u>Native sequence</u>	M227 of the recombinant protein is equivalent to M1 of human DUSP22
<u>Accession number</u>	GenBank NM_020185.

Recombinant DUSP22 amino acid sequence:

```

1  MSPILGYWKI  KGLVQPTRL  LEYLEEKYEE  HLYERDEGDK  WRNKKFELGL  EFPNLPYYID
61  GDVKLTQSMA  IIRYIADKHN  MLGGCPKERA  EISMLEGAVL  DIRYGVSRIA  YSKDFETLKV
121  DFLSKLPEML  KMFEDRLCHK  TYLNGDHVTH  PDFMLYDALD  VVLYMDPMCL  DAFPKLVCFK
181  KRIEAIQID  KYLKSSKYIA  WPLQGWQATF  GGGDHPPKSD  LVPRGSMGNG  MNKILPGLYI
241  GNFKDARDAE  QLSKNKVTHI  LSVHDSARPM  LEGVKYLCIP  AADSPSQNLT  RHFKESIKFI
301  HECRLRGESC  LVHCLAGVSR  SVTLVIAYIM  TVTDFGWEDA  LHTVRAGRSC  ANPNVGFQRQ
361  LQEFEKHEVH  QYRQWLKEEY  GESPLQDAEE  AKNILAAPGI  LKFWAFLRRL

```

Recombinant DUSP22 nucleotide sequence:

```

1  atgtccccta  tactaggtta  ttggaaaatt  aagggccttg  tgcaaccac  tcgacttctt
61  ttggaatata  ttgaagaaaa  atatgaagag  catttgtatg  agcgcgatga  aggtgataaa
121  tggcgaaaca  aaaagtttga  attgggttgg  gagtttccca  atcttcctta  ttatattgat
181  ggtgatgtta  aattaacaca  gtctatggcc  atcatacgtt  atatactga  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  acatggaccc  aatgtgcctg  gatgcgttcc  caaaattagt  ttgttttaaa
541  aaacgtattg  aagctatccc  acaaattgat  aagtacttga  aatccagcaa  gtatatagca
601  tggcctttgc  agggctggca  agccacgttt  ggtggtggcg  accatcctcc  aaaatcggat
661  ctggttccgc  gtggatccat  ggggaatggg  atgaacaaga  tcctgcccgg  cctgtacatc
721  ggcaacttca  aagatgccag  agacgcggaa  caattgagca  agaacaaggt  gacacatatt
781  ctgtctgtcc  acgatagtgc  caggcctatg  ttggaggagg  ttaaataacct  gtgcatccca
841  gcagcggatt  caccatctca  aaacctgaca  agacatttca  aagaaagat  taaattcatt
901  cacgagtgcc  ggctccgctg  tgagagctgc  cttgtacact  gcctggccgg  ggtctccagg
961  agcgtgacac  tggatgatcg  atacatcatg  accgtcactg  actttggctg  ggaggatgcc
1021  ctgcacaccg  tgcgtgctgg  gagatcctgt  gccaaaccca  acgtgggctt  ccagagacag
1081  ctccaggagt  ttgagaagca  tgaggtccat  cagtatcggc  agtggctgaa  ggaagaatat
1141  ggagagagcc  ctttgaggga  tgcagaagaa  gccaaaaaca  ttctggccgc  tccaggaatt
1201  ctgaagttct  gggcctttct  cagaagactg  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.