

## Certificate of Analysis

### Met (D1246H), active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-915, 14-915-K, 14-915M

Parent Lot # D7SN029N

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, Met amino acids 974–end, containing the mutation D1246H, expressed by baculovirus in Sf21 insect cells. Purified using Ni<sup>2+</sup>/NTA agarose.

Met D1246H is a somatic mutation that has been identified in patients with papillary renal carcinomas (HPRC). *In vitro* studies have demonstrated enhanced kinase activity from this mutant. (Schmidt *et al*, (1997), Nature Genetics, **16**: 68-73 and Jeffers *et al* (1997), PNAS, **94**: 11445-11450).

Purity 87% by SDS-PAGE and Coomassie blue staining. MW = 50.1kDa.

**Formulation:** 2.069mg/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.

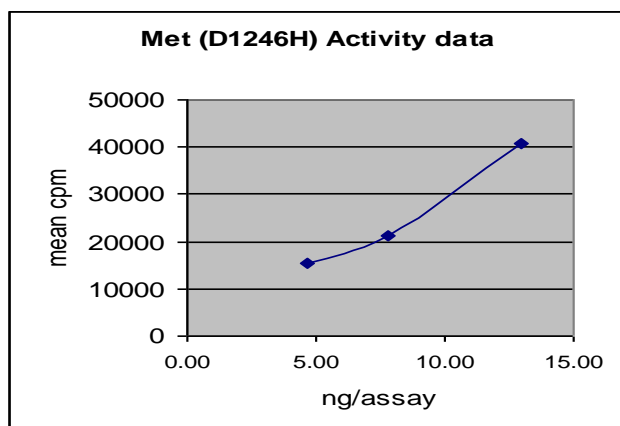
**Specific Activity (Parent lot # D7SN029N):** 1450U/mg, where one unit of Met (D1246H), active activity is defined as 1nmol phosphate incorporated into 250µM (KKKGQEEEEYVFIE) per minute at 30°C with a final ATP concentration of 100µM.

**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.

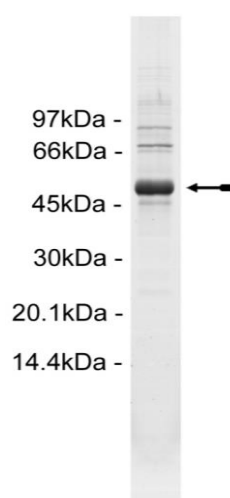
**FOR IN VITRO RESEARCH USE ONLY  
NOT FOR USE IN HUMANS OR ANIMALS**

### Quality Control Testing

**Kinase Assay:** 4.7–12.9ng of this lot of enzyme phosphorylated 250µM (KKKGQEEEEYVFIE) 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 Met 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 Met (D1246H), active.

## Certificate of Analysis

### Kinase Assay Protocol

#### Stock Solutions:

1. **5 x Reaction Buffer:** 40mM MOPS/NaOH pH7.0, 1mM EDTA.
2. **Na<sub>3</sub>VO<sub>4</sub>:** Use at a final assay concentration of 1mM. Prepare a 100mM stock and add 0.25µl of stock per assay point.
3. **Na-β-glycerophosphate:** Use at a final assay concentration of 5mM. Prepare a 1M stock and add 0.125µl of stock per assay point.
4. **(KKKGQEEYVFIE):** Use at a final assay concentration of 250µM. Prepare a 2.5mM stock and add 2.5µl of stock per assay point.
5. **Met (D1246H), active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 4.7–12.9ng per assay point.
6. **[γ-<sup>33</sup>P]ATP:** 2.5 x MgAc/[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 **(KKKGQEEYVFIE)**.
3. Add **2.5µl (4.7–12.9ng) Met (D1246H), active**.
4. Add 4.625µl of dH<sub>2</sub>O.
5. Add 0.25µl Na<sub>3</sub>VO<sub>4</sub>.
6. Add 0.125µl Na-β-glycerophosphate.
7. Add 10µl of diluted [γ-<sup>33</sup>P]ATP mixture.
8. Incubate for 10 minutes at 30°C.
9. Stop the reaction by adding 5µl of 3% phosphoric acid.
10. Transfer a 10µl aliquot onto the appropriate area of a **P30 Filtermat**.
11. Wash the filtermat three times for 5 minutes with 75mM phosphoric acid.
12. Wash the filtermat once for 2 minutes with methanol.
13. Transfer the filtermat to a sealable plastic bag and add 4ml of scintillation cocktail.
14. 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

### Met (D1246H) Sequence Information

<b><u>Protein</u></b>	Human Met
<b><u>Tags</u></b>	N-terminal 6His
<b><u>Native sequence</u></b>	K8 of the recombinant protein is equivalent to K974 of human Met
<b><u>Accession number</u></b>	GenBank J02958. The recombinant protein contains the amino acid substitutions A1209G and V1290L with reference to GenBank J02958. Both substitutions are reported in GenBank BU595386, BQ315895 and BQ316491.

#### **Recombinant Met (D1246H) amino acid sequence:**

```

1 MHHHHHKKR KQIKDLGSEL VRYDARVHTP HLDRLVSARS VSPTTEMVSN ESDYRATFP
61 EDQFPNSSQN GSCRQVQYPL TDMSPILTSQ DSDISSPLLQ NTVHIDLAL NPELVQAVQH
121 VVIGPSSLIV HFNEVIGRGH FGCVYHGTL DNDGKKIHCA VKSLNRITDI GEVSQFLTEG
181 IIMKDFSHPN VLSLLGICLR SEGSPVVLP YMKHGDLRNF IRNETHNPTV KDLIGFGLQV
241 AKGMKYLASK KFHVRDLAAR NCMLDEKFTV KVADFGLARH MYDKEYYSVH NKTGAKLPVK
301 WMALESQTQ KFTTKSDVWS FGVLLWELMT RGAPPYDPVN TFDITVYLLQ GRRLLQPEYC
361 PDPLYEVMLK CWHPKAEMRP SFSELVSRI S AIFSTFIGEH YVHVNATYVN VKCVAPYPSL
421 LSEEDNADDE VDTRPASFE TS
  
```

#### **Recombinant Met (D1246H) nucleotide sequence:**

```

1 atgcatcacc atcaccatca taaaaagcga aagcaaatta aagatctggg cagtgaatta
61 gttcgctacg atgcaagagt acacactcct catttggata ggcttgtaag tgcccgaagt
121 gtaagcccaa ctacagaaat ggtttcaaat gaatctgtag actaccgagc tacttttcca
181 gaagatcagt ttcctaattc atctcagaac ggttcatgcc gacaagtgca gtatcctctg
241 acagacatgt cccccatcct aactagtggg gactctgata tatccagtcc attactgcaa
301 aatactgtcc acattgacct cagtgtctta aatccagagc tgggtccaggc agtgcagcat
361 gtagtgattg ggcccagtag cctgattgtg catttcaatg aagtcatagg aagagggcat
421 tttggttggt tatatcatgg gactttgttg gacaatgatg gcaagaaaat tcaactgtgct
481 gtgaaatcct tgaacagaat cactgcataa ggagaagttt cccaatttct gaccgagga
541 atcatcatga aagattttag tcatcccaat gtcctctcgc tcctgggaat ctgcctgcca
601 agtgaagggt ctccgctggt ggtcctacca tacatgaaac atggagatct tcgaaatttc
661 attcgaaatg agactcataa tccaactgta aaagatccta ttggctttgg tcttcaagta
721 gccaaaggca tgaatatcct tgcaagcaaa aagtttgtcc acagagactt ggctgcaaga
781 aactgtatgc tggatgaaaa attcacagtc aaggttgctg attttggctc tgccagacac
841 atgtatgata aagaatacta tagtgtcac aacaaaacag gtgcaaagct gccagtgaag
901 tggatggctt tggaaagtct gcaaactcaa aagtttacc acaagtcaga tgtgtggtcc
961 tttggcgtgc tcctctggga gctgatgaca agaggagccc caccttatcc tgacgtaaac
1021 accttgata taactgttta cttgttgcaa gggagaagac tcctacaacc cgaatactgc
1081 ccagaccctt tatatgaagt aatgctaaaa tgctggcacc ctaaagccga aatgcgccc
1141 tccttttctg aactggtgtc ccgatatca gcgatcttct ctactttcat tggggagcac
1201 tatgtccatg tgaacgttac ttatgtgaac gtaaaatgtg tcgctccgta tccttctctg
1261 ttgtcatcag aagataacgc tgatgatgag gtggacacac gaccagcctc cttctgggg
1321 acatcatag
  
```

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.