

## Certificate of Analysis

### Met, active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-526, 14-526-K, 14-526M

Parent Lot # WAB0316

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 residues 974–end, expressed by baculovirus in Sf21 insect cells. Purified using Ni<sup>2+</sup>/NTA-agarose. Purity 61% by SDS-PAGE and Coomassie blue staining. MW = 50kDa.

**Specific Activity (Parent lot# WAB0316):** 2683U/mg, where one unit of Met, active activity is defined as 1nmol phosphate incorporated into 250µM (KKKSPGEYVNIEFG) per minute at 30°C with a final ATP concentration of 100µM.

**Formulation:** 0.85mg/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.

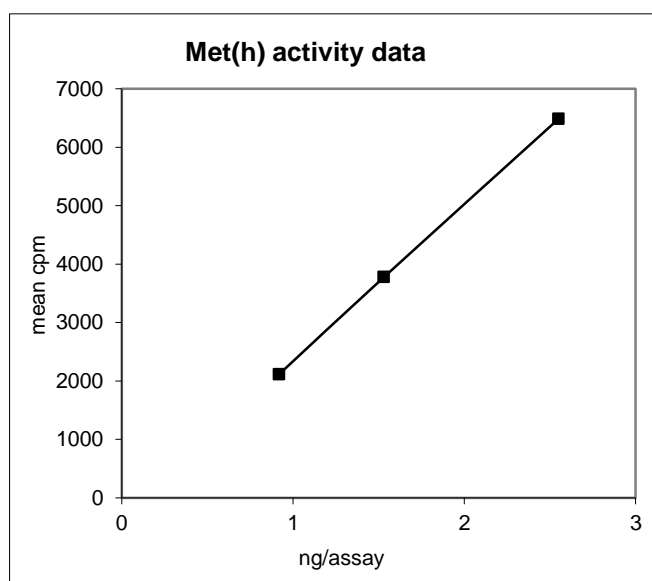
**Storage and Stability:** On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

**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.

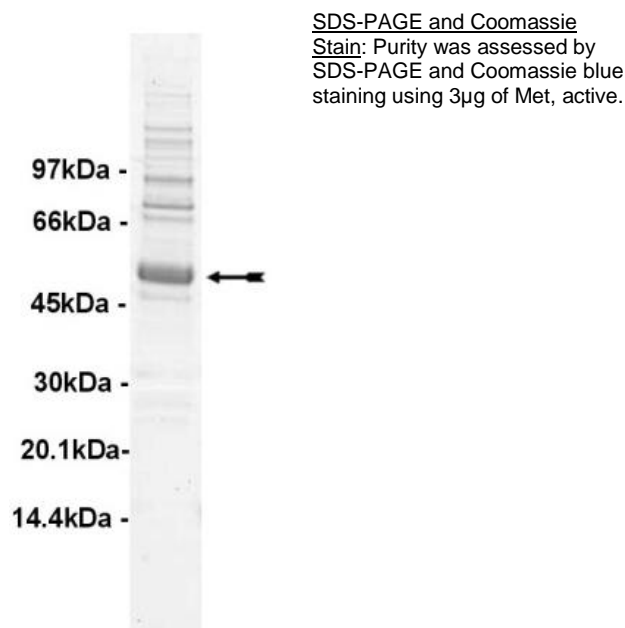
**FOR IN VITRO RESEARCH USE ONLY  
NOT FOR USE IN HUMANS OR ANIMALS**

### Quality Control Testing

**Kinase Assay:** 0.9–2.6ng of this lot of enzyme phosphorylated 250µM in the assay described on page two. Assay background was subtracted from the actual counts to yield the results.



**MS Tryptic Fingerprint:** Confirmed identity Met with the translated 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. **(KKKSPGEYVNIIEFG):** Use at a final assay concentration of 25 $\mu$ M. Prepare a 2.5mM stock and add 2.5 $\mu$ l of stock per assay point.
3. **Met, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 0.9–2.6ng per assay point.
4. **[ $\gamma$ -<sup>33</sup>P]ATP:** 2.5 x magnesium acetate/[ $\gamma$ -<sup>33</sup>P]ATP cocktail: 25mM MgAc and 0.25mM ATP to which is added [ $\gamma$ -<sup>33</sup>P]ATP (specific activity approximately 500 - 800cpm/pmol as required.)

#### Assay Procedure (96 well plate format):

1. Add 5 $\mu$ l of 5 x reaction buffer per assay to wells.
2. Add 2.5 $\mu$ l of **(KKKSPGEYVNIIEFG)**.
3. Add **2.5 $\mu$ l (0.9–2.6ng) Met, active**.
4. Add 5 $\mu$ l of dH<sub>2</sub>O.
5. Add 10 $\mu$ l of diluted [ $\gamma$ -<sup>33</sup>P]ATP mixture.
6. Incubate for 10 minutes at 30°C.
7. Stop the reaction by adding 5 $\mu$ l of 3% phosphoric acid.
8. Transfer a 10 $\mu$ 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 $\mu$ l of 30% phosphoric acid.

## Certificate of Analysis

### Met 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. This construct contains the conflicts A1209G and V1290L with respect to GenBank J02958. Both conflicts are reported in GenBank BU595386, BQ315895 and BQ316491.

#### **Recombinant Met amino acid sequence:**

```

1 MHHHHHHKKR KQIKDLGSEL VRYDARVHTP HLDRLVSARS VSPTTEMVSN ESVDYRATFP
61 EDQFPNSSQN GSCRQVQYPL TDMSPILTSG DSDISSPLLQ NTVHIDL SAL NPELVQAVQH
121 VVIGPSSLIV HFNEVIGRGH FGCVYHG TLL DNDGKKIHCA VKSLNRITDI GEVSQFLTEG
181 IIMKDFSHPN VLSLLGICLR SEGSP LVVLP YMKHGD LRFN IRNETHNPTV KDLIGFGLQV
241 AKGMKYLASK KFVHRDLAAR NCMLDEKFTV KVADFGLARD MYDKEYYSVH NKTGAKLPVK
301 WMALES LQTQ KFTTKSDVWS FGVLLWELMT RGAPPY PDVN TFDITVYLLQ GRRLLQPEYC
361 PDPLYEVMLK CWHPKAEMRP SFSELVSRIS AIFSTFIGEH YVHV NATYVN VKCVAPYPSL
421 LSSEDNADDE VDTRPAS FWE TS
  
```

#### **Recombinant Met 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 cagtgtctca aatccagagc tgggtccaggc agtgcagcat
361 gtagtgattg ggcccagtag cctgattgtg catttcaatg aagtcatagg aagagggcat
421 tttggttggt tatatcatgg gactttgttg gacaatgatg gcaagaaaat tcaactgtgct
481 gtgaaatcct tgaacagaat cactgacata ggagaagttt cccaatttct gaccgaggga
541 atcatcatga aagattttag tcatcccaat gtcctctcgc tcctgggaat ctgcttgcca
601 agtgaagggt ctccgctggt ggtcctacca tacatgaaac atggagatct tcgaaatttc
661 attcgaatg agactcataa tccaactgta aaagatccta ttggctttgg tcttcaagta
721 gccaaaggca tgaatatatc tgcaagcaaa aagtttgctc acagagactt ggctgcaaga
781 aactgtatgc tggatgaaaa attcacagtc aaggttgctg attttggctc tgccagagac
841 atgtatgata aagaatacta tagtgtacac aacaaaacag gtgcaaagct gccagtgaag
901 tggatggcct tggaaagtct gcaaaactca aagtttacc acaagtcaga tgtgtggtcc
961 tttggcgtgc tcctctggga gctgatgaca agaggagccc caccttatcc tgacgtaaac
1021 acctttgata taactgttta cttgttgcaa gggagaagac tcctacaacc cgaatactgc
1081 ccagaccctt tataatgaag aatgctaaaa tgctggcacc ctaaagccga aatgcgccc
1141 tccttttctg aactggtgtc ccggatatca gcgatcttct ctactttcat tggggagcac
1201 tatgtccatg tgaacgctac ttatgtgaac gtaaaatgtg tcgctccgta tccttctctg
1261 ttgtcatcag aagataacgc tgatgatgag gtggacacac gaccagcctc cttctgggag
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.