

## Certificate of Analysis

### Src, active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-326, 14-326-K, 14-326M

Parent Lot # WAB0003

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, full-length, human Src expressed by baculovirus in Sf21 insect cells. Purified using Ni<sup>2+</sup>/NTA-agarose. Purity 95.3% by SDS-PAGE and Coomassie blue staining. MW = 61.7kDa.

**Specific Activity (Parent lot# WAB0003):** 1022U/mg, where one unit of Src activity is defined as 1nmol phosphate incorporated into 250µM cdc2 substrate peptide (cat# 12-140) per minute at 30°C with a final ATP concentration of 100µM.

**Formulation:** 1.128mg/ml of enzyme in 50mM Tris/HCl pH7.5, 150mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 5% glycerol, 1mM benzamidine, 0.2mM PMSF, 0.1% 2-mercaptoethanol. Frozen solution.

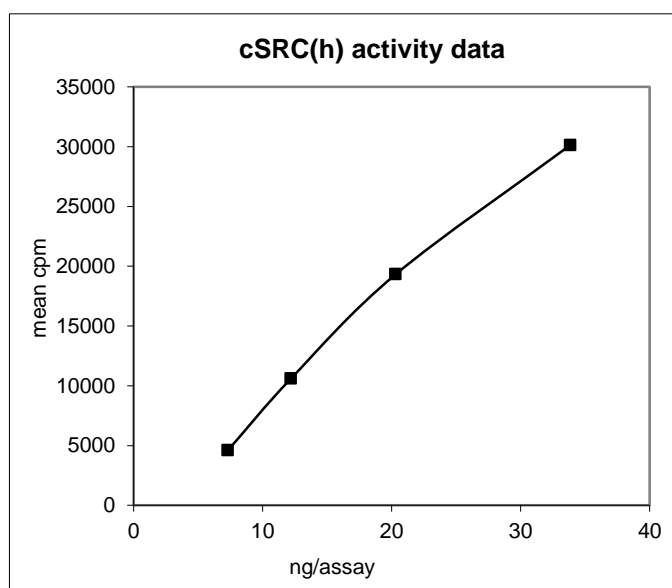
**Storage and Stability:** On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

**Handling Recommendations:** Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled microcentrifuge tubes and immediately snap-freeze the vials in liquid nitrogen prior to re-storage at -70°C.

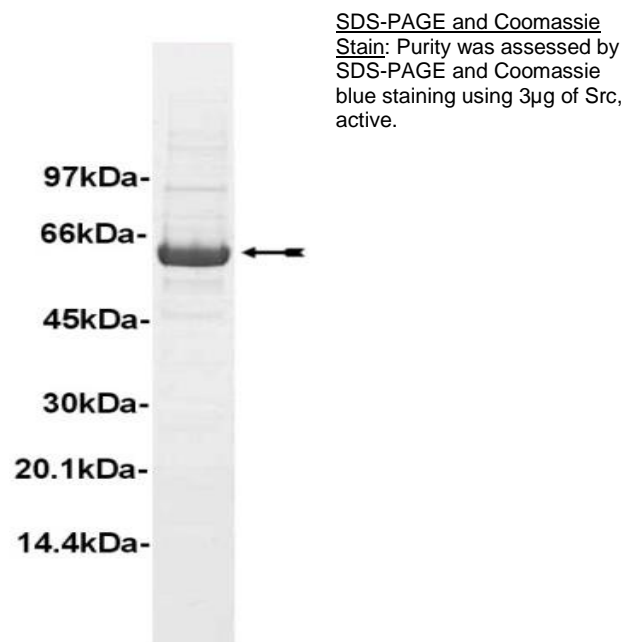
**FOR IN VITRO RESEARCH USE ONLY  
NOT FOR USE IN HUMANS OR ANIMALS**

### Quality Control Testing

**Kinase Assay:** 7–34ng of this lot of enzyme phosphorylated 250µM cdc2 substrate peptide 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 Src with the translated native 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. cdc2 peptide (KVEKIGEGTYGVVYK):** Use at a final assay concentration of 250µM. Prepare a 2.5mM stock and add 2.5µl of stock per assay point.
- 3. Src, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 5% glycerol, 0.01% Brij-35, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 7–34ng 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 5µl of 5 x reaction buffer per assay to appropriate wells.
2. Add 2.5µl of **cdc2 substrate peptide**.
3. Add **2.5µl (7–34ng) Src, active**.
4. Add 5µl of dH<sub>2</sub>O.
5. Add 10µl of diluted [ $\gamma$ -<sup>33</sup>P]ATP mixture.
6. Incubate for 10 minutes at 30°C.
7. Stop the reaction by adding 5µl of 3% phosphoric acid.
8. Transfer a 10µl 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 1µl of 30% phosphoric acid.

## Certificate of Analysis

### Src Sequence Information

<b>Protein</b>	human Src
<b>Tags</b>	N-Terminal 6His
<b>Native sequence</b>	M14 of the recombinant protein is equivalent to M1 of human Src
<b>Accession number</b>	EMBL K03218

#### Recombinant Src amino acid sequence:

```

1 MSYYHHHHH DEFMSNKS PKDASQRRRS LEPAENVHGA GGGAFPASQT PSKPASADGH
61 RGPAAAFAPA AAEPKLFGGF NSSDVTSPQ RAGPLAGGVT TFVALYDYES RTETDLSEFK
121 GERLQIVNNT EGDWWLAHSL STGQTGYIPS NYVAPSDSIQ AEEWYFGKIT RRESERLLL
181 AENPRGTFVL RESETTKGAY CLSVSDFDNA KGLNVKHYKI RKLDSGGFYI TSRTQFNSLQ
241 QLVAYYSKHA DGLCHRLTTV CPTSKPQTQG LAKDAWEIPR ESLRLEVKLG QGCFGEVWVG
301 TWNGTTRVAI KTLKPGTMSP EAFLQEAQVM KCLRHEKLVQ LYAVVSEEPY YIVTEYMSKG
361 SLLDFLKGET GKYLRLPQLV DMAAQIASGM AYVERMNYVH RDLRAANILV GENLVCKVAD
421 FGLARLIEDN EYTARQGAKE PIKWTAPEAA LYGRFTIKSD VWSFGILLTE LTTKGRVPYP
481 GMVNREVLQD VERGYRMPCP PECPELHDL MCQCWRKEPE ERPTFEYLQA FLEDYFTSTE
541 PQYQPGENL
  
```

#### Recombinant Src nucleotide sequence:

```

1 atgtcgtact accatcacca tcaccatcac gatgaattca tgggtagcaa caagagcaag
61 cccaaggatg ccagccagcg gcgccgagc ctggagcccg ccgagaacgt gcacggcgct
121 ggcgggggcg cttccccgc ctcgcagacc cccagcaagc cagcctcggc cgacggccac
181 gcgggcccca gcgggcctt cgccccgcg gccgccgagc ccaagctgtt cggaggcttc
241 aactcctcgg acaccgtcac ctccccgag agggcgggccc cgctggccgg tggagtgacc
301 acctttgtgg ccctctatga ctatgagctc aggacggaga cagacctgtc cttcaagaaa
361 ggcgagcggc tccagattgt caacaacaca gagggagact ggtggctggc ccactcgtc
421 agcacaggac agacaggcta catcccagc aactacgtgg cgccctccga ctccatccag
481 gctgaggagt ggtatcttgg caagatcacc agacgggagt cagagcgggt actgctcaat
541 gcagagaacc cgagagggac cttcctcgtg cgagaaagtg agaccacgaa aggtgcctac
601 tgccctctcag tgtctgactt cgacaacgcc aagggcctca acgtgaagca ctacaagatc
661 cgcaagctgg acagcggcgg cttctacatc acctcccga cccagttcaa cagcctgcag
721 cagctggtgg cctactactc caaacacgcc gatggcctgt gccaccgcct caccaccgtg
781 tgccccacgt ccaagccgca gactcagggc ctggccaagg atgcctggga gatccctcgg
841 gagtcgctgc ggtggagggt caagctgggc cagggtgctc ttggcgagggt gtggatgggg
901 acctggaacg gtaccaccag ggtggccatc aaaaccctga agcctggcac gatgtctcca
961 gaggccttcc tgcaggaggc ccaggatcat aagaagtga ggcatgagaa gctggtgcag
1021 ttgtatgctg tggtttcaga ggagcccatt tacatcgtca cggagtacat gagcaagggg
1081 agtttgctgg actttctcaa gggggagaca ggcaagtacc tgcggctgcc ctacgtggtg
1141 gacatggctg ctcagatcgc ctcaggcatg gcgtacgtgg agcggatgaa ctacgtccac
1201 cgggaccttc gcgagccaa catcctggtg ggagagaacc tgggtgtgcaa agtgccgac
1261 tttgggctgg ctcggctcat tgaagacaat gagtacacgg cgcggcaagg tgccaaattc
1321 cccatcaagt ggacggctcc agaagctgcc ctctatggcc gcttcacat caagtcggac
1381 gtgtggtcct tcgggatcct gctgactgag ctcaccacaa agggacgggt gccctaccct
1441 gggatggtga accgcagggt gctggaccag gtggagcggg gctaccggat gccctgcccg
1501 ccggagtgtc ccgagtcctt gcacgacctc atgtgccagt gctggcggaa ggagcctgag
1561 gagcggccca cttcgagta cctgcaagcc ttctggagg actacttcac gtccaccgag
1621 cccagctacc agcccgggga gaacctctag
  
```

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.