

Certificate of Analysis

Lck, active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-442, 14-442-K, 14-442M

Parent Lot # WAE0125

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 Lck expressed in Sf21 insect cells. Purified using Ni²⁺/NTA agarose. Purity 96% by SDS-PAGE and Coomassie blue staining. MW = 59kDa.

Specific Activity (Parent lot# WAE0125): 394U/mg, where one unit of Lck, active activity is defined as 1nmol phosphate incorporated into 250µM cdc 2 peptide per minute at 30°C with a final ATP concentration of 100µM.

Formulation: 0.99mg/ml of enzyme in 50mM Tris/HCl, pH7.5, 0.1mM EGTA, 150mM NaCl, 1mM benzamidine, 0.03% Brij-35, 0.2mM PMSF, 0.1% 2-mercaptoethanol, 5% glycerol. Frozen solution.

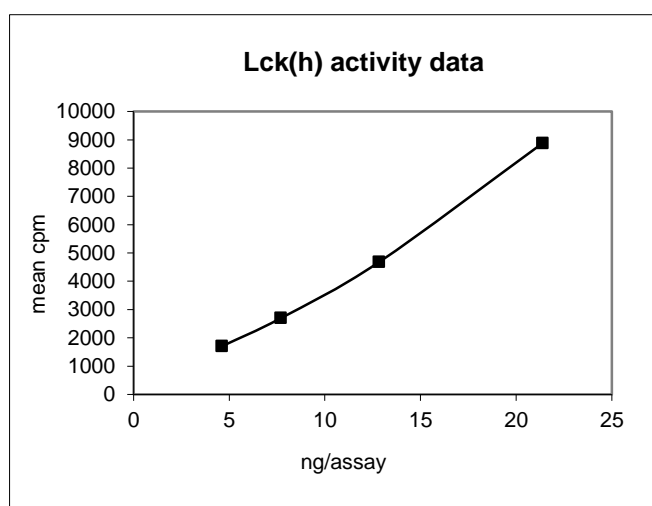
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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.

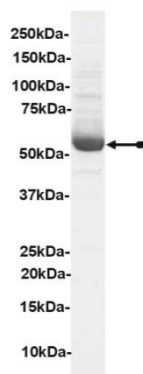
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 4.62–21.38ng of this lot of enzyme phosphorylated 250µM cdc2 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 Lck 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 Lck, active.

Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

1. **4 x Reaction Buffer:** 200mM Tris/HCl pH7.5, 0.4mM EGTA, 0.4mM Na₃VO₄.
2. **cdc2 Substrate Peptide:** Use a final concentration of 250µM. Make a 2.5mM stock. Add 2.5µl of stock per assay point.
3. **Lck, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 5% glycerol, 0.01% Brij-35, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 4.62–21.38ng per assay point.
4. **[γ-³³P]ATP:** 2.5 x magnesium acetate/[γ-³³P]ATP cocktail: 25mM MgAc and 0.25mM ATP to which is added [γ-³³P]ATP (specific activity approximately 500 - 800cpm/pmol as required.)

Assay Procedure (96 well plate format):

1. Add 6.25µl of 4 x reaction buffer per assay to wells.
2. Add 2.5µl of substrate peptide **cdc2**.
3. Add **2.5µl (4.62–21.38ng) Lck, active**.
4. Add 3.75µl of dH₂O.
5. Add 10µl of diluted [γ-³³P]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 twice 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 (background cpm.)

Certificate of Analysis

Lck Sequence Information

Protein	Human Lck
Tags	N-terminal 6His
Native sequence	G9 of the recombinant protein is equivalent to G1 of human Lck
Accession number	Swiss Prot P06239

Recombinant Lck amino acid sequence:

```

1 MHHHHHMGK GCSHPEDDW MENIDVCENC HYPIVPLDGK GTLLIRNGSE VRDPLVTYEG
61 SNPPASPLQD NLVIALHSYE PSHDGLGFE KGEQLRILEQ SGEWKKAQSL TTGQEGFIPF
121 NFVAKANSLE PEPWFFKNLS RKDAERQLLA PGNTHGSFLI RESESTAGSF SLSVRDFDQN
181 QGEVVKHYKI RNLNDGGFYI SPRITFPGLH ELVRHYTNAS DGLCTRLSRP CQTQKPQKPW
241 WEDEWEVPRE TLKLVERLGA GQFGEVVMGY YNGHTKVAVK SLKQGSMSPD AFLAEANLMK
301 QLQHQRLVRL YAVVTQEPIY IITEYMENGS LVDFLKTSPG IKLTINKLLD MAAQIAEGMA
361 FIEERNYIHR DLRAANILVS DTLSCKIADF GLARLIEDNE YTAREGAKFP IKWTAPEAIN
421 YGFTIKSDV WSGILLTEI VTHGRIPYPG MTNPEVIQNL ERGYRMVRPD NCPPELYQLM
481 RLCWKERPED RPTFDYLRV LEDFFTATEG QYQPQP
  
```

Recombinant Lck nucleotide sequence:

```

1 atgcatcacc atcaccatca tatgggctgt ggctgcagct cacaccgga agatgactgg
61 atgaaaaaca tcgatgtgtg tgagaactgc cattatccca tagtcccact ggatggcaag
121 ggacagctgc tcatccgaaa tggctctgag gtgcgggacc cactggttac ctacgaaggc
181 tccaatccgc cggctcccc actgcaagac aacctgggta tcgctctgca cagctatgag
241 ccctctcacg acggagatct gggctttgag aagggggaac agctccgcat cctggagcag
301 agcggcgagt ggtggaaggc gcagtccttg accacgggccc aggaaggctt catccccctt
361 aatthttgtgg ccaaagcgaa cagtctggag cccgaacctt ggttcttcaa gaacctgagc
421 cgcaaggacg cggagcggca gctcctggcg cccgggaaca ctcacggctc cttcctcatc
481 cgggagagcg agagcaccgc gggatcgttt tcaactgtcg tccgggactt cgaccagaac
541 cagggagagg tggtgaaaca ttacaagatc cgtaatctgg acaacgggtg cttctacatc
601 tcccctcgaa tcaactttcc cggcctgcat gaactgggtc gccattacac caatgcttca
661 gatgggctgt gcacacgggt gagccgccc tgccagacc agaagcccc gaagccgtgg
721 tgggaggacg agtgggaggt tcccaggag acgctgaagc tgggtggagc gctgggggct
781 ggacagttcg gggaggtgtg gatgggttac tacaacgggc acacgaaggt ggcggtgag
841 agcctgaagc aggcagcat gtccccggac gccttctgg tccaggccaa cctcatgag
901 cagctgcaac accagcggct ggttcggctc tacgctgtgg tcaccagga cccatctac
961 atcatcactg aatacatgga gaatgggagt ctagtggatt ttctcaagac ccttcaggc
1021 atcaagtga ccatcaaaa actcctggac atggcagccc aaattgcaga aggcattggc
1081 ttattgaag agcgaatta tattcatcgt gacctcggg ctgccaacat tctggtgtct
1141 gacaccctga gctgcaagat tgcagacttt ggcctagcac gcctcattga ggacaacgag
1201 tacacagcca gggagggggc caagtttccc attaagtgga cagcgccaga agccattaac
1261 tacgggacat tcaccatcaa gtcagatgtg tggctctttg ggatcctgct gacggaaatt
1321 gtcaccacag gccgcatccc ttaccagggt atgaccaacc cggaggtgat tcagaacctg
1381 gagcaggct accgcatggt ggccttgac aactgtccag aggagctgta ccaactcatg
1441 aggctgtgct ggaaggagcg cccagaggac cggcccacct ttgactacct gcgcagtgtg
1501 ctggaggact tcttcacggc cacagagggc cagtaccagc ctcagccttg a
  
```

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.