

Certificate of Analysis

Abl (mouse), active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-459, 14-459-K, 14-459M

Parent Lot # LB164P86

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 mouse Abl, amino acids 27–end, expressed by baculovirus in Sf21 insect cells. Purified using Ni²⁺/NTA agarose. Purity 61% by SDS-PAGE and Coomassie blue staining. MW = 123.5kDa.

Specific Activity (Parent lot# LB164P86): 1195U/mg, where one unit of Abl (mouse), active activity is defined as 1nmol phosphate incorporated into 50µM Abltide (EAIYAAPFAKKK) per minute at 30°C with a final ATP concentration of 100µM.

Formulation: 3.19mg/ml of enzyme in 50mM Tris/HCl pH7.5, 150mM NaCl, 0.1mM EGTA, 0.03% Brij-35, 270mM sucrose, 0.2mM PMSF, 1mM benzamidine, 0.1 % 2-mercaptoethanol. Frozen solution.

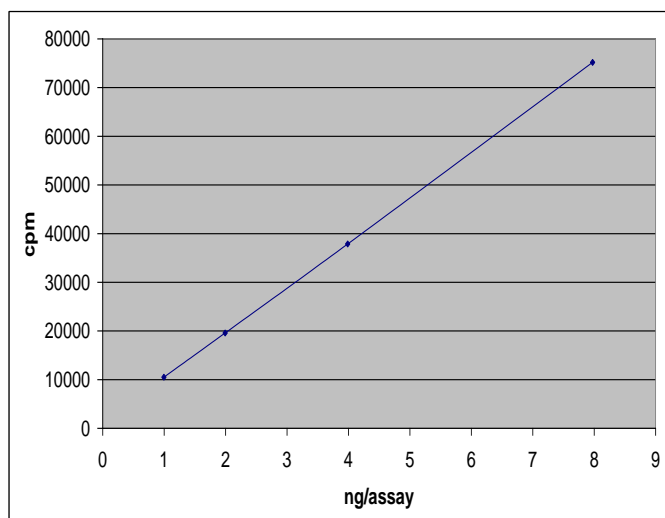
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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: 1–8ng of this lot of enzyme phosphorylated 50µM Abltide 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 Abl (mouse) with the translated sequence listed on pages three and four.

97 kDa

66 kDa

45 kDa

30 kDa

20.1 kDa

kDa

4.4 kDa



SDS-PAGE and Coomassie Stain: Purity was assessed by SDS-PAGE and Coomassie blue staining using 3µg of Abl (mouse), active.

Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

1. **5 x Reaction Buffer:** 40mM MOPS/NaOH pH7.0, 1mM EDTA.
2. **Abltide (EAIYAAPFAKKK):** Use at a final assay concentration of 50 μ M. Prepare a 500 μ M stock and add 2.5 μ l of stock per assay point.
3. **Abl (mouse), active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 1–8ng 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 5 μ l of 5 x reaction buffer per assay to wells.
2. Add 2.5 μ l of **Abltide (EAIYAAPFAKKK)**.
3. Add **2.5 μ l (1–8ng) Abl (mouse), active**.
4. Add 5 μ l of dH₂O.
5. Add 10 μ l of diluted [γ -³³P] ATP mixture.
6. Incubate for 10 minutes at 30°C.
7. Stop the reaction by adding 5 μ l of 3% phosphoric acid.
8. Transfer a 10 μ 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 μ l of 30% phosphoric acid.

Certificate of Analysis

Mouse Abl Sequence Information

<u>Protein</u>	mouse Abl
<u>Tags</u>	N-Terminal 6His
<u>Native sequence</u>	E30 of the recombinant protein is equivalent to E27 of mouse Abl
<u>Accession number</u>	GenBank J02995. The recombinant protein contains four amino acids which conflict with J02995, these are L782P, G784R, W785L, and L786V, and are reported in GenBank XM_130089.

Recombinant Mouse Abl amino acid sequence:

```

1 MSYYHHHHHH DYDIPTTENL YFQGAMDPGE ALQRPVASDF EPQGLSEEAR WNSKENLLAG
61 PSENDPNLHV ALYDFVASGD NTLSITKGEK LRVLGYNHNG EWCEAQTKNG QGWVPSNYIT
121 PVNSLEKHSW YHGPFVSRNAA EYLLSSGING SFLVRESESS PGQRSISLRY EGRVYHYRIN
181 TASDGKLYVS SESRFNTLAE LVHHHSTVAD GLITTLHYPA PKRKNKPTIYG VSPNYDKWEM
241 ERTDITMKHK LGGGQYGEVY EGVWKKYSLT VAVKTLKEDT MEVEEFLKEA AVMKKEIKHPN
301 LVQLLGVCTR EPPFYIITEF MTYGNLLDYL RECNRQEVSA VVLLYMATQI SSAMEYLEKK
361 NFIHRDLAAR NCLVGENHLV KVADFGLSRL MTGDTYTAHA GAKFPIKWTA PESLAYNKFS
421 IKSDVWAFGV LLWEIATYGM SPYPGIDLSQ VYELLEKDYR MERPEGCPEK VYELMRACWQ
481 WNPSDRPSFA EIHQAFETMF QESSISDEVA KELGKRGRTRG GAGSMLQAPE LPTKTRTCRR
541 AAEQKDAPDT PELLHTKGLG ESDALDSEPA VSPLLPRKER GPPDGSLNED ERLLPDRDKT
601 NLFSAIILKK KKMAPTTPKR SSSFREMDDQ PDRRGASEDD SRELGNPPA LTSDAAEPTK
661 SPKASNGAGV PNGAFREPGN SGFRSPHMWK KSSTLTGSRL AAAEEESGMS SSKRFLRSCS
721 ASCMPHGARD TEWRSVTLPR DLPSAGKQFD SSTFGGHKSE KPALPRKRTS ESRSEQVAKS
781 TAMPPPRLVK KNEEAAEEGF KDTESSPGSS PPSLTPKLLR RQVTASPSSG LSHKEEATKG
841 SASGMGTPAT AEPAPPSNKV GLSKASSEEM RVRRHKHSSE SPGRDKGRLA KLKPAAPPPP
901 ACTGKAGKPA QSPSQEAGEA GGPTTKKCTS LAMDAVNTPD TKAGPPGEGE RKPVPSPVVK
961 PQSTAKPPGT PTSPVSTPST APAPSPLAGD QQPSSAAFIP LISTRVSLRK TRQPPERIAS
1021 GTITKGVVLD STEALCLAIS RNSEQMASHS AVLEAGKNLY TFCVSYVDSI QQMRNKFAFR
1081 EAINKLESNL RELQICPATA SSGPAATQDF SKLLSSVKEI SDIVRR

```

Recombinant Mouse Abl nucleotide sequence:

```

1 atgtcgtact accatcacca tcaccatcac gattacgata tcccaacgac cgaaaacctg
61 tattttcagg gcgccatgga tccggggcgaa gccctgcaga ggccagtggc atctgacttt
121 gagccccagg gtctcagcga agcagctcga tggaaactcca agggaaacct tcttgctggg
181 cccagtgaat atgaccccaa cttttttgtg gcactctatg attttgtggc cagtggagat
241 aacactctca gcatcactaa aggtgaaaag ctccgggtct tgggttataa tcacaatggg
301 gaatggtgtg aagcccaaac gaaaaatggc caagatggg tcccaagcaa ctacatcacc
361 cccgtcaaca gctgggagaa acattctctg tatcatggcc ctgtatctcg gaatgctgct
421 gagtatctgc tgagcagcgg aatcaacggc agcttcttag tgcgggagag tgagagtagc
481 cctggccaga gatccatctc gctgcggtat gaagggaggg tgtaccacta caggatcaac
541 actgcctctg atggcaagct gtacgtgtcc tccgagagcc gcttcaaac tctggctgag
601 ttagttcacc atcactccac ggtggctgat ggcctcatca ccacactcca ctaccagct
661 cccaagcgca acaagcccac tatctacggt gtgtcccca actacgaca gtgggaaatg
721 gagcgcaccg acatcaccat gaagcacaag ttgggtggag gccagtagcg ggaggtgtac
781 gagggcggtt ggaagaagta cagcctcact gtggcctgta agacctgaa ggaggacacc
841 atggaggtgg aggagttcct gaaggaagcg gcggtgatga aggagatcaa acaccctaac
901 ctggtgcagc tgctaggggt gtgtaccgga gaaccacat tctacataat cactgagttc
961 atgacctatg ggaacctgct ggactacctg agggagtgtg accggcagga ggtgagcgcc
1021 gtggtactgc tctacatggc cacacagatc tcatcagcca tggagtactt ggagaagaag
1081 aacttcatcc acagagacct tgctgcccgg aactgctgg taggggaaaa cacttgggtg
1141 aaggtggctg attttggcct gagcaggttg atgacagggg acacctacac ggcccatgct
1201 ggagccaaat tccccatcaa atggaccgca cctgagagcc tggcctacaa caagttctcc

```

Certificate of Analysis

```

1261 atcaagtcgg acgtgtgggc atttgagta ttgctctggg agattgctac ctatggcatg
1321 tcaccttacc cgggaattga cctgtctcag gtttatgagc tgctggaaaa agactaccgc
1381 atggagcgcc ctgaaggctg cccggagaag gtctacgagc tcatgcgagc atgttggcag
1441 tggaaaccct ctgaccggcc ctcccttctg gaaatccacc aagcctttga aacctatgtc
1501 caggaatcca gtatctcaga tgaggctggg aaggagctgg ggaaacgagg cacgagagga
1561 ggtgctggga gtatgctgca ggcccagag ctgcccacca agaccagaac ctgcaggaga
1621 gcagctgagc agaaagatgc gcctgacacc cctgagctgc tccacacgaa gggcctggga
1681 gaaagcgatg cactggacag tgagcctgct gtatcgccac tgcttcctcg gaaagagcgc
1741 gggcccccag acggcagcct aaatgaagat gagcgccttc tccccagaga cagaaagacc
1801 aacctgttca gcgctttgat caagaagaag aagaaaatgg cgccgacgcc ccctaagcgc
1861 agcagttcct tccgagagat ggatggccag ccagaccgca gaggggctag tgaggatgac
1921 agcagggaac tctgcaatgg accaccagct ctcacctcag acgcagcaga gcctaccaag
1981 tccccaaagg ccagcaatgg ggctggcgct cctaattggag ccttcgggga gccgggcaac
2041 tcaggcttcc gttctcccca catgtggaaa aagtccagca cactgaccgg gagccgcctg
2101 gctgctgccg aagaggagag cggcatgagc tccagtaagc gcttcctgcg ttcttgctcg
2161 gcctcctgca tgccccatgg ggcaagggac acagagtggc ggtcggctac gctgcctcga
2221 gacctgccgt ctgctggcaa gcagtttgac tcatccacct ttggagggca caaaagcgaa
2281 aagccagctc tgccctcggaa acgcaccagt gagagcaggt ctgagcaggt ggccaaaagc
2341 acggcgatgc cccctccccg gctggtgaag aagaacgagg aggctgctga agaaggcttc
2401 aaagacacag aatccagccc tggctccagc cctcccagct tgactcccaa actcctccgc
2461 aggcaggtca ctgcctctcc ttctcttggc ctctctcaca aggaagaggc caccaagggc
2521 agtgctcag gcgatggggac tccggccact gcagagccag cccccccag caacaaagtg
2581 ggctcagca aggcctcctc tgaggagatg cgcgtaagga ggcacaagca cagctcggag
2641 tccccagggg gagacaaggg gcgactggct aagctcaagc ctgccccgcc gcctcctcct
2701 gcctgcacag gaaaagcagg caagcccgca cagagcccca gccaaagagg cggggaggga
2761 ggggggcccc caaagacaaa atgcacgagt ctggctatgg atgctgtgaa cactgacccc
2821 accaaggccc gccacactgg agaaggactg agaaagcctg tgccccatc tgtgccaaag
2881 ccccagtcga cggctaagcc tccagggact cccaccagcc cggctctccac cccctccaca
2941 gcaccagctc cttcaccctt ggctggggac cagcagccat cttctgcccgc cttcatcccc
3001 ctcatatcaa cccgtgtgtc tcttaggaag acccgccagc cgccagagcg cattgcccagt
3061 ggcaccatca ccaaggggtg ggttctggac agtactgagg ccctgtgctt tgccatctcc
3121 cggaactcag agcagatggc cagccacagt gctgtactgg aggctggcaa gaacctgtac
3181 actttctgtg tgagctatgt ggactctatc cagcagatga ggaacaagtt tgccttccgt
3241 gaggctatca acaagctgga gagcaacctc cgagagctgc agatctgccc tgccacagcc
3301 tccagtgggc cagctgccac ccaagacttc agcaagctgc tcagctctgt gaaggagatc
3361 agcgacattg tccggaggta g

```

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.