

Certificate of Analysis

Abl (M351T), active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-757, 14-757-K, 14-757M

Parent Lot # 1643914

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 Abl, residues 27–end containing the M351T mutation. Expressed by baculovirus in Sf21 insect cells. Purified using Ni²⁺/NTA agarose.

The Abl tyrosine kinase inhibitor STI571 is an effective therapy for stable-phase chronic myeloid leukaemia patients. Many patients responding to STI571 later relapse due to a reactivation of Bcr-Abl activity. In certain cases this appears to correlate with the presence of the M351T mutation, which confers resistance to drug binding (Shah et al., (2002), Cancer Cell 2, 117-125).

Purity 40% by SDS-PAGE and Coomassie staining. MW = 121.4kDa.

Formulation: 0.783mg/ml of enzyme in 50mM Tris/HCl pH7.5, 150mM NaCl, 270mM sucrose, 1mM benzamidine, 0.2mM PMSF, 0.1mM EGTA, 0.1% 2-mercaptoethanol, 0.03% Brij-35. 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.

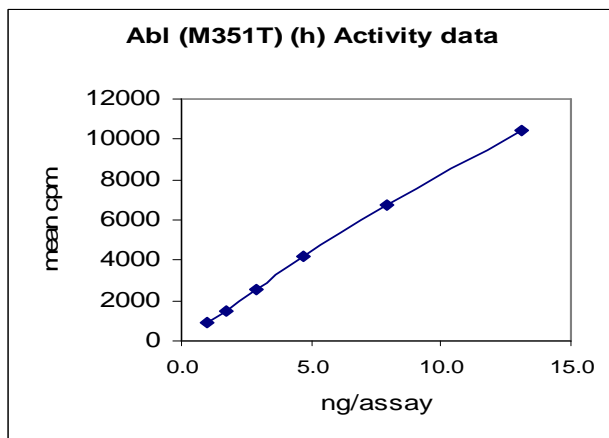
Specific Activity (Parent lot# 1643914): 899U/mg, where one unit of Abl (M351T) activity is defined as 1nmol phosphate incorporated into 50µM Abltide (EAIYAAPFAKCK) per minute at 30°C with a final ATP concentration of 100µM.

Handling Recommendations: Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled 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: 1.0–13.1ng of this lot of enzyme phosphorylated 50µM Abltide (EAIYAAPFAKCK) 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 product identity as Abl (M351T) 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. **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 (M351T), active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 1.0–13.1ng 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**.
3. Add **2.5 μ l (1.0–13.1ng Abl (M351T), 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

Abl (M351T) Sequence Information

Protein	Human Abl (27–end, M351T)
Tags	N-terminal 6His
Native sequence	E10 of recombinant sequence is equivalent to E27 of native human Abl
Accession number	GenBank U07563

Recombinant Abl (M351T) amino acid sequence:

```

1 MHHHHHHEFE ALQRPVASDF EPQGLSEAAR WNSKENLLAG PSENDPNLFV ALYDFVASGD
61 NTLISITKGEK LRVLGYNHNG EWCEAQTNG QGWVPSNYIT PVNSLEKHSW YHGPVSRNAA
121 EYLLSSGING SFLVRESESS PGQRSISLRY EGRVYHYRIN TASDGKLYVS SESRFNTLAE
181 LVHHHSTVAD GLITTLHYPA PKRKNKPTVYG VSPNYDKWEM ERTDITMKHK LGGGQYGEVY
241 EGVWKKYSLT VAVKTLKEDT MEVEEFLKEA AVMKIKHPN LVQLLGVCTR EPPFYIITEF
301 MTYGNLLDYL RECNRQEVNA VVLLYMATQI SSATEYLEKK NFIHRDLAAR NCLVGENHLV
361 KVADFGLSRL MTGDTYTAHA GAKFPIKWTA PESLAYNKFS IKSDVWAFGV LLWEIATYGM
421 SPYPGIDLSQ VYELLEKDYR MERPEGCPEK VYELMRACWQ WNPSDRPSFA EIHQAFETMF
481 QESSISDEVE KELGKQVVRG AVSTLLQAPE LPTKTRTSRR AAHRDRTDV PEMPHSKGQG
541 ESDPLDHEPA VSPLLPRKER GPPEGGLNED ERLLPKDKKT NLFSALIKKK KKTAPTTPKR
601 SSSFREMDGQ PERRGAGEEE GRDISNGALA FTPLDTADPA KSPKPSNGAG VPNGALRESG
661 GSGFRSPHLW KKSSTLTSSR LATGEEEGGG SSSKRFLRSC SASCVPHGAK DTEWRSVTLF
721 RDLQSTGRQF DSSTFGGHKS EKPALPRKRA GENRSDQVTR GTVTPPRRLV KKNEEADEV
781 FKDIMESSPG SSPPNLTPKP LRRQVTVAPA SGLPHKEEAG KGSALGTPAA AEPVTPTSKA
841 GSGAPGGTSK GPAEESRVRH HKHSSESPGR DKGKLSRLKP APPPPPAASA GKAGGKPSQS
901 PSQEAAGEAV LGAKTKATSL VDAVNSDAK PSQPGGLKK PVLPATPKPQ SAKPSGTPIS
961 PAPVPSTLPS ASSALAGDQP SSTAFAIPLIS TRVSLRKRTRQ PPERIASGAI TKGVVLDSTE
1021 ALCLAISRNS EQMASHSAVL EAGKNLYTFC VSYVDSIQQM RNKFVAFREAI NKLENNLREL
1081 QICPATAGSG PAATQDFSKL LSSVKEISDI VQR

```

Recombinant Abl (M351T) nucleotide sequence:

```

1 atgcatcatt accatcacca tgaattcgaa gcccttcagc ggccagtagc atctgacttt
61 gagcctcagg gtctgagtga agccgctcgt tggaaactcca agggaaacct tctcgctgga
121 cccagtgaaa atgaccccaa ccttttcggt gcactgtatg attttgtggc cagtggagat
181 aacactctaa gcataactaa aggtgaaaag ctccgggtct taggctataa tcacaatggg
241 gaatggtgtg aagcccaaac caaaaatggc caaggctggg tccaagcaa ctacatcacg
301 ccagtcaaca gtctggagaa acactcctgg taccatgggc ctgtgtcccg caatgccgct
361 gagtatctgc tgagcagcgg gatcaatggc agcttcttgg tgcgtgagag tgagagcagt
421 cctggccaga ggtccatctc gctgagatac gaaggagggg tgtaccata caggatcaac
481 actgcttctg atggcaagct ctacgtctcc tccgagagcc gcttcaaac cctggccgag
541 ttggttcatt atcattcaac ggtggccgac gggctcatca ccacgctcca ttatccagcc
601 ccaaagcgca acaagcccac tgtctatggt gtgtcccca actacgacaa gtgggagatg
661 gaacgcacgg acatcaccat gaagcacaag ctgggcccgg gccagtagcg ggaggtgtac
721 gagggcgtgt ggaagaaata cagcctgacg gtggccgtga agacctgaa ggaggacacc
781 atggagggtg aagagtctt gaaagaagct gcagtcatga aagagatcaa acaccctaac
841 ctggtgcagc tccttggggc ctgcaccgag gagccccgct tctatatcat cactgagttc
901 atgacctagc ggaacctcct ggactacctg agggagtgca accggcagga ggtgaacgcc
961 gtggtgctgc tgtacatggc cactcagatc tcgtcagcca cggagtacct ggagaagaaa
1021 aacttcatcc acagagatct tgctgcccga aactgcctgg taggggagaa ccacttggtg
1081 aaggtagctg attttggcct gagcaggttg atgacagggg acacctacac agcccatgct
1141 ggagccaagt tccccatcaa atggactgca cccgagagcc tggcctacaa caagttctcc
1201 atcaagtccg acgtctgggc atttggagta ttgcttggg aaattgctac ctatggcatg
1261 tccccttacc cgggaattga cctgtcccag gtgtatgagc tgctagagaa ggactaccgc
1321 atggagcgcc cagaaggctg cccagagaag gtctatgaac tcatgtagc atgttggcag

```

Certificate of Analysis

```

1381  tggaaatccct  ctgaccggcc  ctcctttgct  gaaatccacc  aagcctttga  aacaatgttc
1441  caggaatcca  gtatctcaga  cgaagtggaa  aaggagctgg  ggaaacaagg  cgtccgtggg
1501  gctgtgagta  ccttgctgca  ggccccagag  ctgcccacca  agacgaggac  ctccaggaga
1561  gctgcagagc  acagagacac  cactgacgtg  cctgagatgc  ctactccaa  gggccaggga
1621  gagagcgatc  ctctggacca  tgagcctgcc  gtgtctccat  tgctccctcg  aaaagagcga
1681  ggtcccccg  agggcggcct  gaatgaagat  gagcgccttc  tccccaaaga  caaaaagacc
1741  aacttgttca  gcgccttgat  caagaagaag  aagaagacag  cccaacccc  tccccaaagc
1801  agcagctcct  tccgggagat  ggacggccag  ccggagcgca  gaggggccc  cgaggaagag
1861  ggccgagaca  tcagcaacgg  ggcactggct  ttcaccccct  tggacacagc  tgaccagcc
1921  aagtcccaa  agcccagcaa  tggggctggg  gtcccacatg  gagccctccg  ggagtccggg
1981  ggctcaggct  tccggtctcc  ccacctgtgg  aagaagtcca  gcacgtgac  cagcagccgc
2041  ctagccaccg  gcgaggagga  gggcggtggc  agctccagca  agcgttccct  gcgctcttgc
2101  tccgcctcct  gcgttccca  tggggccaag  gacacggagt  ggaggtcagt  cacgctgcct
2161  cgggacttgc  agtccacggg  aagacagttt  gactcgtcca  catttggagg  gcacaaaagt
2221  gagaagccgg  ctctgcctcg  gaagagggca  ggggagaaca  ggtctgacca  ggtgaccgca
2281  ggcacagtaa  cgcctcccc  caggctgggt  aaaaagaatg  aggaagctgc  tgatgaggtc
2341  ttcaaagaca  tcatggagtc  cagcccgggc  tccagcccgc  ccaacctgac  tccaaaacc
2401  ctccggcggc  aggtcacctg  ggcccctgcc  tcgggcctcc  cccacaagga  agaagctgga
2461  aagggcagtg  ccttagggac  cctgctgca  gctgagccag  tgacccccac  cagcaaagca
2521  ggctcaggtg  caccagggg  caccagcaag  ggcccgcgc  aggagtccag  agtgaggagg
2581  cacaagcact  cctctgagtc  gccagggagg  gacaagggga  aattgtccag  gctcaaacct
2641  gccccgcgc  ccccaccagc  agcctctgca  gggaaaggctg  gaggaaagcc  ctgcagagc
2701  ccgagccagg  aggcggccgg  ggaggcagtc  ctgggcgcaa  agacaaaagc  cagcagctg
2761  gttgatgctg  tgaacagtga  cgctgccaag  cccagccagc  cgggagagg  cctcaaaaag
2821  cccgtgctcc  cggccactcc  aaagccacag  tccgccaagc  cgtcggggag  ccccatcagc
2881  ccagcccccg  ttccctccac  gttgccatca  gcatcctcgg  ccctggcagg  ggaccagccg
2941  tcttccaacc  cttcatccc  totcatatca  acccgagtgt  ctcttcggaa  aaaccgccag
3001  cctccagagc  ggatcgccag  cggcgccatc  accaagggcg  tggctctgga  cagcaccgag
3061  gcgctgtgcc  tcgccatctc  taggaactcc  gagcagatgg  ccagccacag  cgcagtgctg
3121  gaggccggca  aaaacctcta  cacgttctgc  gtgagctatg  tggattccat  ccagcaaagt
3181  aggaacaagt  ttgccttccg  agaggccatc  aacaaactgg  agaataatct  ccgggagctt
3241  cagatctgcc  cggcgacagc  aggcagtggt  ccagcggcca  ctcaggactt  cagcaagctc
3301  ctcagttcgg  tgaaggaaat  cagtgatata  gtgcagaggt  ag

```

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.