

Certificate of Analysis

Abl (Y253F), active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-759, 14-759-K, 14-759M

Parent Lot # 1615549

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 Y253F mutation. Expressed by baculovirus in Sf21 insect cells. Purified using immobilised metal affinity chromatography.

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 Y253F mutation, which confers resistance to drug binding (Shah et al., (2002), Cancer Cell 2, 117-125).

Purity 62% by SDS-PAGE and Coomassie staining. MW = 121.4kDa.

Specific Activity (Parent lot# 1615549): 2351U/mg, where one unit of Abl (Y253F), active activity is defined as 1nmol phosphate incorporated into 50µM Abltide (EAIYAAPFAKCK) per minute at 30°C with a final ATP concentration of 100µM.

Formulation: 1.071mg/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.

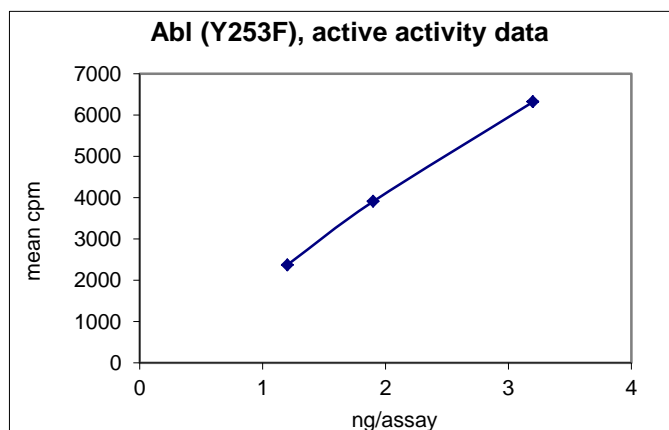
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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.

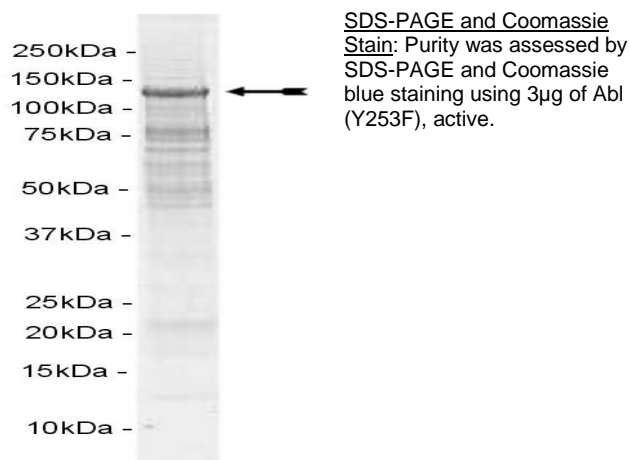
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 1.2–3.2ng 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 (Y253F) 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 (Y253F), active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 0.01% Brij-35, 5% glycerol, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 1.2–3.2ng 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 (1.2–3.2ng) Abl (Y253F), 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 (Y253F) Sequence Information

<u>Protein</u>	Human Abl (27-end, Y253F)
<u>Tags</u>	N-terminal 6His
<u>Native sequence</u>	E10 of recombinant sequence is equivalent to E27 of native human Abl
<u>Accession number</u>	GenBank U07563

Recombinant Abl (Y253F) amino acid sequence:

```

1  MHHHHHEFE  ALQRPVASDF  EPQGLSEEAR  WNSKENLLAG  PSENDPNLFV  ALYDFVASGD
61  NTLSTITKGEK  LRVLGYNHNG  EWCEAQTKNG  QGWVPSNYIT  PVNSLEKHSW  YHGPVSRNAA
121  EYLLSSGING  SFLVRESESS  PGQRSISLRY  EGRVYHYRIN  TASDGKLYVS  SESRFNTLAE
181  LVHHHSTVAD  GLITTLHYPA  PKRNKPTVYG  VSPNYDKWEM  ERTDITMKHK  LGGGQFGEVY
241  EGVWKKYSLT  VAVKTLKEDT  MEVEEFLKEA  AVMKEIKHPN  LVQLLGVCTR  EPPFYIITEF
301  MTYGNLLDYL  RECNRQEVNA  VLLYMATQI  SSAMEYLEKK  NFIHRDLAAR  NCLVGENHLV
361  KVADFGLSRL  MTGDTYTAHA  GAKFPIKWTA  PESLAYNKFS  IKSDVWAFGV  LLWEIATYGM
421  SPYPGIDLSQ  VYELLEKDYR  MERPEGCPEK  VYELMRACWQ  WNPSDRPSFA  EIQAFETMF
481  QESSISDEVE  KELGKQGVRG  AVSTLLQAPE  LPTKTRTSRR  AAEHRDSTDV  PEMPHSKGQG
541  ESDPLDHEPA  VSPLLPRKER  GPPEGGLNED  ERLLPKDKKT  NLFSAIKKK  KKTAPTPPKR
601  SSSFREMDGQ  PERRGAGEEE  GRDISNGALA  FTPLDTADPA  KSPKPSNGAG  VPNGALRESG
661  GSGFRSPHLW  KKSSTLTSSR  LATGEEEGGG  SSKRFLRSC  SASCVPHGAK  DTEWRSVTLF
721  RDLQSTGRQF  DSSTFGGHKS  EKPALPRKRA  GENRSDQVTR  GTVTPPPRLV  KKNEEADEV
781  FKDIMESSPG  SPPNLTPKP  LRRQVTVAPA  SGLPHKEEAG  KGSALGTPAA  AEPVTPSKA
841  GSGAPGGTSK  GPAESRVR  HKHSSSESPGR  DKGKLSRLKP  APPPPAASA  GKAGGKPSQS
901  PSQEAAGEAV  LGAKTKATSL  VDAVNSDAAK  PSQPGEGLKK  PVLPATPKPQ  SAKPSGTPIS
961  PAPVPSTLPS  ASSALAGDQP  SSTAFLPLIS  TRVSLRKTRQ  PPERIASGAI  TKGVVLDSTE
1021  ALCLAISRNS  EQMASHSAVL  EAGKNLYTFC  VSYVDSIQQM  RNKFAFREAI  NKLENNLREL
1081  QICPATAGSG  PAATQDFSKL  LSSVKEISDI  VQR

```

Recombinant Abl (Y253F) nucleotide sequence:

```

1  atgcatcatc  accatcacca  tgaattcgaa  gcccttcagc  ggccagtagc  atctgacttt
61  gagcctcagg  gtctgagtga  agccgctcgt  tggaactcca  aggaaaacct  tctcgtctga
121  cccagtgaaa  atgaccccaa  cttttcgtt  gcactgtatg  attttgtggc  cagtggagat
181  aacactctaa  gcataactaa  aggtgaaaag  ctccgggtct  taggctataa  tcacaatggg
241  gaatggtgtg  aagcccaaac  caaaaatggc  caaggctggg  tccaagcaa  ctacatcacg
301  ccagtcaaca  gtctggagaa  aactcctgg  taccatgggc  ctgtgtcccg  caatgccgct
361  gagtatctgc  tgagcagcgg  gatcaatggc  agcttcttgg  tgcgtgagag  tgagagcagt
421  cctggccaga  ggtccatctc  gctgagatac  gaaggagggg  tgtaccatta  caggatcaac
481  actgcttctg  atggcaagct  ctacgtctcc  tccgagagcc  gcttcaacac  cctggccgag
541  ttggttcatc  atcattcaac  ggtggccgac  gggctcatca  ccacgctcca  ttatccagcc
601  ccaaagcgca  acaagccac  tgtctatggt  gtgtcccca  actacgacaa  gtgggagatg
661  gaagcagcgg  acatcaccat  gaagcacaag  ctgggcgggg  gccagttcgg  ggaggtgtac
721  gagggcgtgt  ggaagaaata  cagcctgacg  gtggccgtga  agacctttaa  ggaggacacc
781  atggaggtgg  aagagttctt  gaaagaagct  gcagtcatga  aagagatcaa  acaccctaac
841  ctggtgcagc  tccttgggg  ctgcaccgg  gagccccgt  tctatatcat  cactgagttc
901  atgacctacg  ggaacctcct  ggactacctg  agggagtgca  accggcagga  ggtgaacgcc
961  gtggtgctgc  tgtacatggc  cactcagatc  tcgtcagcca  tggagtacct  ggagaagaaa
1021  aacttcatcc  acagagatct  tgctgccga  aactgcctgg  taggggagaa  ccacttgggtg
1081  aaggtagctg  attttgcct  gagcaggttg  atgacagggg  acacctacac  agcccatgct
1141  ggagccaagt  tccccatcaa  atggactgca  cccgagagcc  tggcctacaa  caagttctcc
1201  atcaagtccg  acgtctgggc  atttggagta  ttgctttggg  aaattgctac  ctatggcatg

```

Certificate of Analysis

```

1261 tcccccttacc cgggaattga cctgtcccag gtgtatgagc tgctagagaa ggactaccgc
1321 atggagcgc cagaaggctg cccagagaag gtctatgaac tcatgcgagc atgttggcag
1381 tggaatccct 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 ggccttgat caagaagaag aagaagacag cccaacccc tcccaaacgc
1801 agcagctcct tccgggagat ggacggccag ccggagcgca gggggccgg cgaggaagag
1861 ggccgagaca tcagcaacgg ggcactggct ttcaccctc tggacacagc tgaccagcc
1921 aagtccccaa agcccagcaa tggggctggg gtcccaatg gagcctccg ggagtccggg
1981 ggctcaggct ccacgtctcc ccactgtgg aagaagtcca gcacgctgac gcagagccgc
2041 ctagccaccg gcgaggagga tggcggtggc agctccagca agcgcttct cgctcttgc
2101 tccgctcctc gcttccccca tggggccaag gacacggagt ggaggtcagt cacgctgcct
2161 cgggacttgc agtccacggg aagacagttt gactcgtcca catttggagg gcacaaaagt
2221 gagaagccgg ctctgcctcg gaagagggca ggggagaaca ggtctgacca ggtgaccga
2281 ggcacagtaa cgctcccc caggctggtg aaaaagaatg aggaagctgc tgatgaggtc
2341 ttcaaagaca tcatggagtc cagcccgggc tccagcccgc ccaacctgac tccaaaacc
2401 ctccggcggc aggtcacctg ggccccctgc tcgggcctcc cccacaagga agaagctgga
2461 aaggcgagt ccttagggac ccctgctgca gctgagccag tgacccccac cagcaaagca
2521 ggctcaggtg caccaggggg caccagcaag ggccccgccc aggagtccag agtgaggagg
2581 cacaagcact cctctgagtc gccagggagg gacaagggga aattgtccag gctcaaact
2641 gccccgccc ccccaccagc agcctctgca ggggaaggctg gaggaaagcc ctgcagagc
2701 ccgagccagg agggcggcgg ggaggcagtc ctgggcgcaa agacaaaagc cacgagtctg
2761 gttgatgctg tgaacagtga cgctgccaag cccagccagc cgggagaggg cctcaaaaag
2821 cccgtgctcc cggccactcc aaagccacag tccgccaagc cgtcggggac ccccatcagc
2881 ccagccccg ttcctccac gttgccatca gcatcctcgg ccctggcagg ggaccagccg
2941 tcttccaccg ccttcatccc tctatatca acccgagtgt ctcttcggaa aaccgcccag
3001 cctccagagc ggatgccag cggcgccatc accaaggcg tggtcctgga cagcaccgag
3061 gcgctgtgcc tcgccatctc taggaactcc gagcagatgg ccagccacag cgcagtgtg
3121 gaggccggca aaaactcta cacgttctgc gtgagctatg tggattccat ccagcaaatg
3181 aggaacaagt ttgccttccg agaggccatc aacaaactgg agaataatct cggggagctt
3241 cagatctgcc cggcgacagc aggcagtggg 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