

## ABL1 M351T

ABL proto-oncogene 1, non-receptor tyrosine kinase

Recombinant Human Active Protein Kinase

HGNC Symbol: ABL1

Synonyms: c-Abl, JTK7, p150

Product No.: 0990-0000-1

Lot: 006

**Description:** Human ABL1, internal fragment, amino acids P<sub>118</sub>-S<sub>535</sub> (as in [NCBI/Protein](#) entry NP\_005148.2), M<sub>351</sub>T mutant, N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

**Product identity:** ABL1 M351T Lot 006, was confirmed as ABL1 with a mutation M351T by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>Fusion Protein</sub>:** 76,137 Da

**Expression host:** Sf9 insect cells

**Purification:** GST-Affinity Chromatography

**Activation:** This kinase was not activated by special procedures

**Storage buffer:** 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20 % glycerol

**Storage temperature:** -80°C

For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

**Protein concentration:** 0.156 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

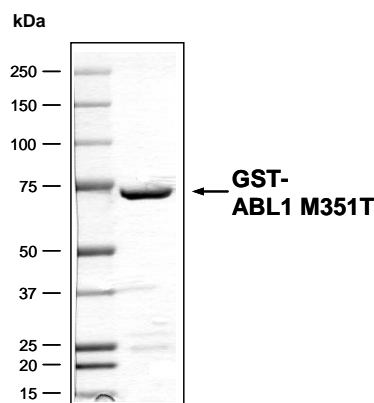
### Biochemical Parameters:

Specific kinase activity (P<sub>i</sub> transfer): 36 pmol/µg × min  
ATP-K<sub>M</sub>: 0.3 µM

### Additional assay technology:

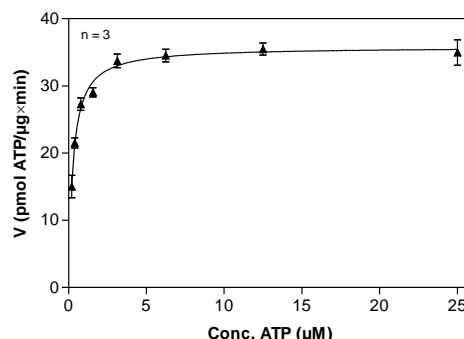
ABL1 M351T Lot 006 was also successfully tested by ProQinase for the use with the ADP-Glo™ Kinase assay from Promega  
ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

### ABL1 M351T Lot 006: Coomassie stain



2.0 µg GST-ABL1 M351T

### ABL1 M351T Lot 006: Determination of V<sub>max</sub> and K<sub>M</sub> value for ATP



- Assay conditions:  
60 mM HEPES-NaOH, pH 7.5  
3 mM MgCl<sub>2</sub>  
3 mM MnCl<sub>2</sub>  
3 µM Na-orthovanadate  
1.2 mM DTT  
50 µg/ml PEG<sub>20,000</sub>  
ATP (variable)  
Substrate: Poly(Ala,Glu;Lys;Tyr)<sub>6:2:5:1</sub>, 20µg/ml  
Kinase: 2 µg/ml
- Filter binding assay  
MSFC membrane (Millipore)

# ABL1 M351T

Product No.: 0990-0000-1

GST-ABL1 M351T Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMG <b>HHHHHG</b>	RDS <b>LEVLFCG</b>	240
241	<b>PLAMLPVNSL</b>	<b>EKHSWYHGPV</b>	<b>SRNAAEYLLS</b>	<b>SGINGSFLVR</b>	<b>ESESSPGQRS</b>	<b>ISLRYEGRVY</b>	300
301	<b>HYRINTASDG</b>	<b>KLYVSSSESRF</b>	<b>NTLAEVLVHHH</b>	<b>STVADGLITT</b>	<b>LHYPAPKRNK</b>	<b>PTVYGVSPNY</b>	360
361	<b>DKWEMERTDI</b>	<b>TMKHKLGGGQ</b>	<b>YGEVYEGVWK</b>	<b>KYSLTVAVKT</b>	<b>LKEDTMEVEE</b>	<b>FLKEAAVMKE</b>	420
421	<b>IKHPNLVQLL</b>	<b>GVCTREPPFY</b>	<b>IITEFMTYGN</b>	<b>LLDYLRECNR</b>	<b>QEVNAVLLY</b>	<b>MATQISSATE</b>	480
481	<b>YLEKKNFIHR</b>	<b>DLAARNCLVG</b>	<b>ENHLVKVADF</b>	<b>GLSRLMTGDT</b>	<b>YTAHAGAKFP</b>	<b>IKWTAPESLA</b>	540
541	<b>YNKFSIKSDV</b>	<b>WAFGVLLWEI</b>	<b>ATYGMSPYPG</b>	<b>IDLSQVYELL</b>	<b>EKDYRMERPE</b>	<b>GCPEKVYELM</b>	600
600	<b>RACWQWNP</b>	<b>RPSFAEIHQA</b>	<b>FETMFQESSI</b>	<b>SDEVEKELGK</b>	<b>QGVRGAVSTL</b>	<b>LQAPELPTKT</b>	660
661	<b>RTS</b>						720

1-218: GST **Red**: HIS6-tag **Green**: 3C cleavage site **blue**: ABL1 fragment **boxed**: M351T

ABL1 wt <sup>1</sup> Amino Acid Sequence							
1	MLEICLKLVG	CKSKKGLSSS	SSCYLEEALQ	RPVASDFEPQ	GLSEAAWNS	KENLLAGPSE	60
61	NDPNLNFVALY	DFVASGDNTL	SITKGEKLRV	LGYNHNGEWC	EAQTKNGQGW	VPSNYIT <b>PVN</b>	120
121	<b>SLEKHSWYHG</b>	<b>PVSRNAAEYL</b>	<b>LSSGINGSFL</b>	<b>VRESESSPGQ</b>	<b>RSISLRYEGR</b>	<b>VYHYRINTAS</b>	180
181	<b>DGKLVSSSE</b>	<b>RFNTLAELVH</b>	<b>HHSTVADGLI</b>	<b>TTLHYPAPKR</b>	<b>NKPTVYGVSP</b>	<b>NYDKWEMERT</b>	240
241	<b>DITMKHKLGG</b>	<b>GQYGEVYEGV</b>	<b>WKKYSLTVAV</b>	<b>KTLKEDTMEV</b>	<b>EEFLKEAAVM</b>	<b>KEIKHPNLVQ</b>	300
301	<b>LLGVCTREPP</b>	<b>FYIITEFMTY</b>	<b>GNLLDYLREC</b>	<b>NRQEVNAVVL</b>	<b>LYMATQISSA</b>	<b>MEYLEKKNFI</b>	360
361	<b>HRDLAARNCL</b>	<b>VGENHLVKVA</b>	<b>DFGLSRLMTG</b>	<b>DTYTAHAGAK</b>	<b>FPIKWTAPES</b>	<b>LAYNKFSIKS</b>	420
421	<b>DVWAFGVLLW</b>	<b>EIATYGMSPY</b>	<b>PGIDLSQVYE</b>	<b>LLEKDYRMER</b>	<b>PEGCPEKVYE</b>	<b>LMRACWQWNP</b>	480
481	<b>SDRPSFAEIH</b>	<b>QAFETMFQES</b>	<b>SISDEVEKEL</b>	<b>GKQGVRGAVS</b>	<b>TLLQAPELPT</b>	<b>KTRTSRRAAE</b>	540
541	HRDITDVPPEM	PHSKGQGESD	PLDHEPAVSP	LLPRKERGPP	EGGLNEDERL	LPKDKKTNLF	600
600	SALIKKKKKT	APTPPKRSSH	FREMDGQPER	RGAGEEEGRD	ISNGALAFTP	LDTADPAKSP	660
661	KPSNGAGVPN	GALRESGGSG	FRSPHLWKKK	STLTSSRLAT	GEEEGGGSSS	KRFLRSCSAS	720
721	CVPHGAKDTE	WRSVTLPRDL	QSTGRQFDSS	TFGGHKSEKP	ALPRKRAGEN	RSDQVTRGTV	780
781	TPPPRLVKKN	EAAAEVFKD	IMESSPGSSP	PNLTPKPLRR	QVTVAPASGL	PHKEEAGKGS	840
841	ALGTPAAAEF	VTPTSKAGSG	APGGTSKGP	EESRVRHKKH	SSESPGRDKG	KLSRLKPAPP	900
901	PPPAASAGKA	GGKPSQSPSQ	EAAGEAVLGA	KTKATSLVDA	VNSDAAKPSQ	PGEGLKPPVL	960
961	PATPKPQSAK	PSGTPISPAP	VPSTLPSASS	ALAGDQPSST	AFIPLISTRV	SLRKTRQPPE	1020
1021	RIASGAIKTK	VVLDSTEALC	LAISRNSEQM	ASHSAVLEAG	KNLYTFCVSY	VDSIQQMRNK	1080
1081	FAFREAINKL	ENNLRELQIC	PATAGSGPAA	TQDFSKLLSS	VKEISDIVQR		1140

**blue**: ABL1 sequence expressed in recombinant protein **Red**: variant in recombinant protein

<sup>1</sup>[NCBI/Protein](https://www.ncbi.nlm.nih.gov/Protein/NP_005148.2) accession number NP\_005148.2

This product was manufactured at ProQinase in Freiburg, Germany, and is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.