

## ProQinase™ ALK G1202R

ALK receptor tyrosine kinase

Recombinant Human Active Protein Kinase

HGNC Symbol: ALK

Synonyms: n/a

Product No.: 1369-0000-1

Lot: 003

**Description:** Human ALK G1202R, internal fragment, amino acids L<sub>1066</sub>-S<sub>1437</sub> (as in [NCBI/Protein](#) entry NP\_004295.2), G1202R mutation, N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

**Product identity:** ALK G1202R Lot 001, was confirmed as ALK by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>Fusion Protein</sub>:** 70,558 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

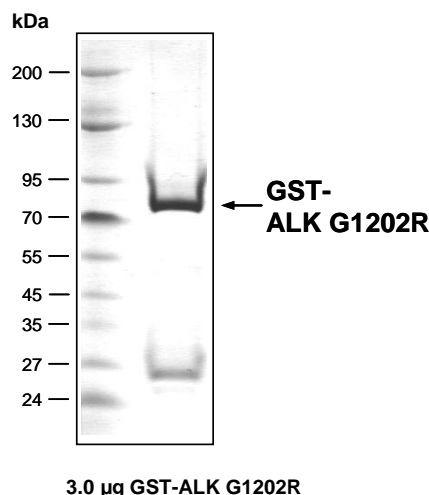
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

**Protein concentration:** 0.108 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

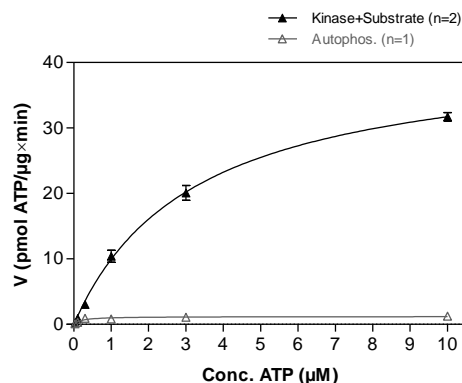
### Biochemical Parameters:

Specific kinase activity (P<sub>i</sub> transfer): 42 pmol/µg × min  
ATP-K<sub>M</sub>: 3.2 µM

### ALK G1202R Lot 003: Coomassie stain



### ALK G1202R Lot 003: 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: TRK-C derived peptide 80 µg/ml  
Kinase: 2 µg/ml
- Filter binding assay  
MSPH membrane (Millipore)

### Additional assay technology:

ALK G1202R Lot 003 was also successfully tested by Reaction Biology 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

## ProQinase™ ALK G1202R

Product No.: 1369-0000-1

GST-ALK G1202R Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMG <b>HHHHHG</b>	RDS <b>LEVLFCG</b>	240
241	<b>PLAMVLQAMQ</b>	<b>MELQSP EYKL</b>	<b>SKLRTSTIMT</b>	<b>DYNPNYCFAG</b>	<b>KTSSISDLKE</b>	<b>VPRKNITLIR</b>	300
301	<b>GLGHGAFGEV</b>	<b>YEQVSGMPN</b>	<b>DPSPLQVAVK</b>	<b>TLPEVCSEQD</b>	<b>ELDFLMEALI</b>	<b>ISKFNHQNIV</b>	360
361	<b>RCIGVSLQSL</b>	<b>PRFILLEMA</b>	<b>GGDLKSFLRE</b>	<b>TRPRPSQPSS</b>	<b>LAML DLLHVA</b>	<b>RDIACGCQYL</b>	420
421	<b>EENHF IHRDI</b>	<b>AARNCLL TCP</b>	<b>GPGRVAKIGD</b>	<b>FGMARDIYRA</b>	<b>SYRKGCCAM</b>	<b>LPVKWMPPEA</b>	480
481	<b>FMEGIFTSKT</b>	<b>DTWSFGVLLW</b>	<b>EIFSLGYMPY</b>	<b>PSKSNQEVLE</b>	<b>FVTSGGRMDP</b>	<b>PKNCPGPVYR</b>	540
541	<b>IMTQCWQHQP</b>	<b>EDRPNFAILL</b>	<b>ERIEYCTQDP</b>	<b>DVINTALPIE</b>	<b>YGPLVEEEEK</b>	<b>VPVRPKDPEG</b>	600
601	<b>VPPLLV SQQA</b>	<b>KREEERS</b>					660

1-218: GST **Red**: HIS6-tag **Green**: 3C cleavage site **blue**: ALK fragment **boxed**: G1202R

ALK wt <sup>1</sup> Amino Acid Sequence							
1	MGAIGLLWLL	PLLSTAAVG	SGMGTGQRAG	SPAAGPPLQP	REPLSYSRLQ	RKSLAVDFV	60
61	PSLFRVYARD	LLLPPSSSEL	KAGRPEARGS	LALDCAPLLR	LLGPAPGVSW	TAGSPAPAEA	120
121	RTLSRVLKG	SVRKLRRAKQ	LVLELGEEAI	LEGCVGPPGE	AAVGLLQFNL	SELSFWWIRQ	180
181	GEGRLRIRLM	PEKKASEVGR	EGRLSAAIRA	SQPRLLFQIF	GTGHSSLESP	TNMPPSPDY	240
241	FTWNLTWIMK	DSFPFLSHRS	RYGLECSFDF	PCELEYSPL	HDLRNQSWSW	RRIPSEEASQ	300
301	MDLLDGP GAE	RSKEMPRGSF	LLLNTSADSK	HTILSPWMS	SSEHCTLA VS	VHRHLQPSGR	360
361	YIAQLLP HNE	AAREILLMPT	PGKHGWTVLQ	GRIGRPD NPF	RVALEYISSG	NRSLSAVDF	420
421	ALKNCSEGTS	PGSKMALQSS	FTCWNGTVLQ	LGQACDFHQD	CAQGEDESQM	CRKLPVGFYC	480
481	NFEDGFCGWT	QGTLSPTHPQ	WQVRTLK DAR	FQDHQDHALL	LSTTDVPASE	SATVTSATFP	540
541	APIKSSPCEL	RMSWLIRGVL	RGNVSLV LVE	NKTGKEQGRM	VWHVAAYEGL	SLWQWMV LPL	600
600	LDVSDRFWLQ	MVAWWGQGSR	AIVAFDNISI	SLDCYLTISG	EDKILQNTAP	KSRNLFERNP	660
661	NKELKPGENS	PRQTPIFDPT	VHWFLLTTCGA	SGPHGPTQAQ	CNNAYQNSNL	SVEVGSEGPL	720
721	KGIQIWKVPA	TDTYSISGYG	AAGGKGKNT	MMRSHGVS VL	GIFNLEKDDM	LYILVGQQGE	780
781	DACPSTNQLI	QKVCIGENNV	I EEEIRVNRS	VHEWAGGGGG	GGGATYVFKM	KDGVVPLII	840
841	AAGGGGRAYG	AKTDTFHPER	LENNSSVLGL	NGNSGAAGGG	GGWNDNTSLL	WAGKSLQEGA	900
901	TGGHSCPQAM	KKWGWETRGG	FGGGGGGCSS	GGGGGGYIGG	NAASNNDPEM	DGEDGVSFIS	960
961	PLGILYTPAL	KVMEGHGEVN	IKHYLNCSHC	EVDECHMDPE	SHKVICFCDH	GTVLAEDGVS	1020
1021	CIVSPTPEPH	LPLSLILSVV	TSALVAALVL	AFSGIMIVYR	RKHQELQAMQ	<b>MELQSP EYKL</b>	1080
1081	<b>SKLRTSTIMT</b>	<b>DYNPNYCFAG</b>	<b>KTSSISDLKE</b>	<b>VPRKNITLIR</b>	<b>GLGHGAFGEV</b>	<b>YEQVSGMPN</b>	1140
1141	<b>DPSPLQVAVK</b>	<b>TLPEVCSEQD</b>	<b>ELDFLMEALI</b>	<b>ISKFNHQNIV</b>	<b>RCIGVSLQSL</b>	<b>PRFILLEMA</b>	1200
1201	<b>GGDLKSFLRE</b>	<b>TRPRPSQPSS</b>	<b>LAML DLLHVA</b>	<b>RDIACGCQYL</b>	<b>EENHF IHRDI</b>	<b>AARNCLL TCP</b>	1260
1261	<b>GPGRVAKIGD</b>	<b>FGMARDIYRA</b>	<b>SYRKGCCAM</b>	<b>LPVKWMPPEA</b>	<b>FMEGIFTSKT</b>	<b>DTWSFGVLLW</b>	1320
1321	<b>EIFSLGYMPY</b>	<b>PSKSNQEVLE</b>	<b>FVTSGGRMDP</b>	<b>PKNCPGPVYR</b>	<b>IMTQCWQHQP</b>	<b>EDRPNFAILL</b>	1380
1381	<b>ERIEYCTQDP</b>	<b>DVINTALPIE</b>	<b>YGPLVEEEEK</b>	<b>VPVRPKDPEG</b>	<b>VPPLLV SQQA</b>	<b>KREEERS</b>	1440
1441	PPPLPTTSSG	KAACKPTAAE	ISVRVPRGPA	VEGGHV NMAF	SQSNPPSELH	KVHGSRNKPT	1500
1501	SLWNPTYGSW	FTEKPTKKN	PIAKKEPHDR	GNLGLGESC T	VPPNVATGRL	PGASLLLEPS	1560
1561	SLTANMKEVP	LFRLRHFP CG	NVNYGYQQQG	LPLEAATAPG	AGHYEDTILK	SKNSMNQPGP	1620

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

<sup>1</sup>[NCBI/Protein](#) accession number NP\_004295.2