

ProQinase™ NPM1ALK

Nucleophosmin anaplastic lymphoma kinase fusionprotein

Recombinant Human Active Protein Kinase

HGNC Symbol: n/a

Synonyms: n/a

Product No.: 1236-0000-1

Lot: 005

Description: Human pathological fusionprotein NPM1ALK, full length, amino acids M₁-P₆₈₀ (as in NCBI/Protein entry AAA58698.1), untagged, expressed in Sf9 insect cells

Product identity: NPM1ALK Lot 005, was confirmed as NPM1ALK by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 75,695 Da

Expression host: Sf9 insect cells

Expression: Baculovirus infected Sf9 cells

Purification: GST-Affinity Chromatography, followed by 3C mediated removal of the GST tag

Activation: This kinase was not activated by special procedures

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

Storage temperature: -80°C

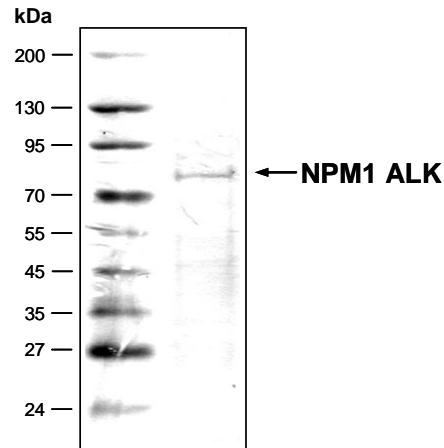
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.189 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:
Specific kinase activity (P_i transfer): 145 pmol/µg×min
ATP-K_M: 5.2 µM

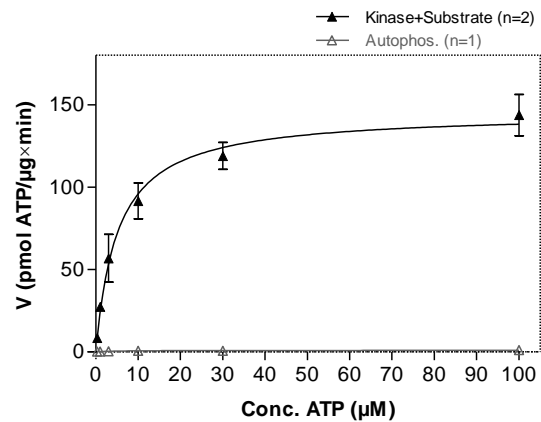
Additional assay technology: NPM1ALK Lot 005 was also successfully tested by Reaction Biology for the use with the ADP-Glo™ Kinase assay from ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

NPM1ALK Lot 005: Coomassie stain



2.0 µg NPM1 ALK

NPM1ALK Lot 005: Determination of V_{max} and K_M value for ATP



Determination of K_M value & Specific activity:

- Assay conditions:
 - 60 mM HEPES-NaOH, pH 7.5
 - 3 mM MgCl₂
 - 3 mM MnCl₂
 - 3 µM Na-orthovanadate
 - 1.2 mM DTT
 - 50 µg / ml PEG_{20,000}
 - ATP (variable)
 - Substrate: TRK-C derived peptide
 - NPM1 ALK: 1.0 µg / ml
- Filter binding assay
 - MSPH membrane (Millipore)

Recombinant Proteins



ProQinase™ NPM1ALK

Product No.: 1236-0000-1

NPM1ALK Recombinant Fusion Protein Amino Acid Sequence								
1	GPLA	MEDSMD	MDMSPLRPQN	YLFGCELKAD	KDYHFKVDND	ENEHQLSLRT	VSLGAGAKDE	60
61	LHIVEAEAMN	YEGSPIKVTL	ATLKMSVQPT	VSLGGFEITP	PVVLRLKCGS	GPVHISGQHL		120
121	VVYRRKHQEL	QAMQMELOSP	EYKLSKLRTS	TIMTDYNPNY	CFAGKTSSIS	DLKEVPRKNI		180
181	TLIRGLGHGA	FGEVYEGQVS	GMPNDPSPLQ	VAVKTLPEVC	SEQDELDFLM	EALIISKFNH		240
241	QNIVRCIGVS	LQSLPRFILL	ELMAGDLKS	FLRETRPRPS	QPSSLAMLDL	LHVARDIACG		300
301	CQYLEENHFI	HRDIAARNCL	LTCPGPGRVA	KIGDFGMARD	IYRASYYRKG	GCAMPLVKWM		360
361	PPEAFMEGIF	TSKTDTWSFG	VLLWEIFSLG	YMPYPSKSNQ	EVLEFVTSGG	RMDPPKNCPG		420
421	PVYRIMTQCW	QHQPEDRPNF	AIILERIEYC	TQDPDVINTA	LPIEYGPLVE	EEEKVPVRPK		480
481	DPEGVPLLIV	SQQAQKREER	SPAAPPPLPT	TSSGKAAKPK	TAAEVSVRVP	RGPAVEGGHV		540
541	NMAFSQSNPP	SELHIVHGSR	NKPTSLWNPT	YGSWFTEKPT	KKNNPIAKKE	PHERGNGGLE		600
601	GSCVPPNVA	TGRLPGASLL	LEPSSLTANM	KEVPLFRLRH	FPCGNVNYGY	QQQGLPLEAA		660
661	TAPGAGHYED	TILKSKNSMN	QPGP					720

1-4: legacy of 3C cleavage blue: NPM1ALK boxed: variation from RefSeq

NPM1ALK wt ¹ amino acid sequence							
1	MEDSMDMDMS	PLRPQNYLFG	CELKADKDYH	FKVDNDENEH	QLSLRTVSLG	AGAKDELHIV	60
61	EAEAMNYEGS	PIKVTLATLK	MSVQPTVSLG	GFEITPPVVL	RLKCGSGPVH	ISGQHLVVYR	120
121	RKHQELQAMQ	MELQSPEYKL	SKLRSTIMT	DYNPNYCFAG	KTSSISDLKE	VPRKNITLIR	180
181	GLGHGAFGEV	YEGQVSGMPN	DPSPLQVAVK	TLPEVCSEQD	ELDFLMEALI	ISKFNHQNIV	240
241	RCIGVSLQSL	PRFILLELMA	GGDLKSFLRE	TRPRPSQPSS	LAMLDDLHVA	RDIACGCQYL	300
301	EENHFIHRDI	AARNCLLTCP	GPGRVAKIGD	FGMARDIYRA	SYRKGGCAM	LPVKWMPPEA	360
361	FMEGIFTSKT	DTWSFGVLLW	EIFSLGYMPY	PSKSNQEVLE	FVTSGGRMDP	PKNCPGPVYR	420
421	IMTQCWQHQP	EDRPNFMAIL	ERIEYCTQDP	DVINTALPIE	YGPLVEEEEK	VPVRPKDPEG	480
481	VPLLIVSQQA	KREERSPAA	PPPLPTTSSG	KAARKPTAAE	VSVRVPRGPA	VEGGHVNMAF	540
541	SQSNPPSELH	KVHGSRNKPT	SLWNPTYGSW	FTEKPTKNN	PIAKKEPHDR	GNLGLEGSCT	600
601	VPPNVATGRL	PGASLLLEPS	SLTANMKEVP	LFRLRHFPCC	NVNYGYQQQG	LPLEAATAPG	660
661	AGHYEDTILK	SKNSMNQPGP					720

bold letters: expressed part of NPM1 (blue) and ALK (green) RED letters: variant in Fusionprotein

¹NCBI/Protein accession number AAA58698.1

K551R and D589E: SNP variations see NCBI/dbSNP IDs: rs1881420, rs1881421