

ProQinase™ EML4 ALK F1174L

Echinoderm microtubule-associated protein-like 4 anaplastic lymphoma kinase fusionprotein

Recombinant Human Active Protein Kinase

HGNC Symbol: n/a

Synonyms: n/a

Product No.: 1247-0000-1

Lot: 003

Description: Human pathological fusionprotein EML4 ALK, full length, amino acids M₁-P₁₀₅₉ (as in NCBI/Protein entry BAF73611.1) with a F1174L mutation, N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: EML4 ALK F1174L Lot 003, was confirmed as EML4 ALK by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 145.413 Da

Expression: Baculovirus infected Sf9 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
Avoid repeated freeze-thaw cycles!

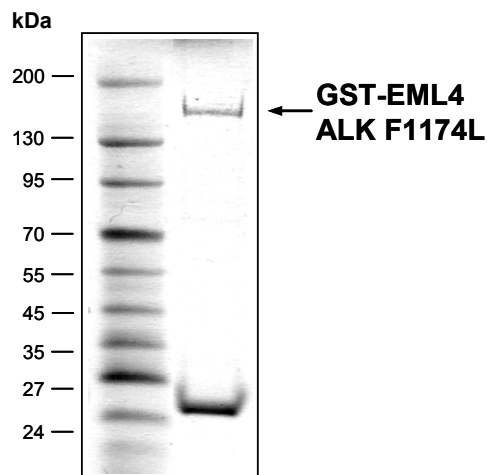
Protein concentration: 0.111 µg/µl (Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

Specific activity: 29 pmol/µg×min

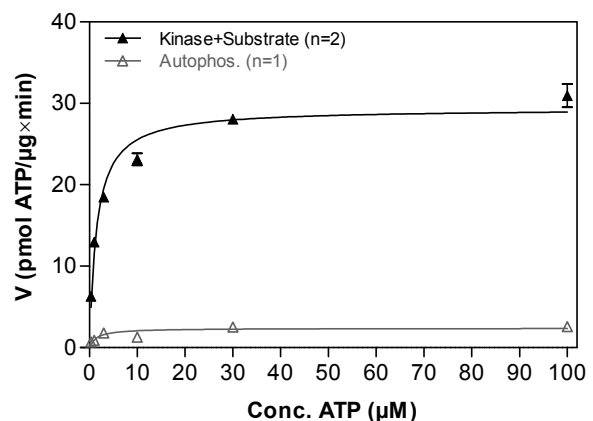
ATP-K_M: 1.5 µM

**EML4 ALK F1174L Lot 003:
Coomassie stain**



2.0 µg GST-EML4 ALK F1174L

**EML4 ALK F1174L Lot 003:
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
 - EML4 ALK F1174L: 2.0 µg / ml
- Filter binding assay
- MSPH membrane (Millipore)

Additional assay technology: EML4 ALK F1174L Lot 003

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



Recombinant Proteins

ProQinase™ EML4 ALK F1174L

Product No.: 1247-0000-1

Recombinant Proteins

EML4 ALK F1174L Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQ SMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQG WQATF	GGGDHPPKSD	PMGHHHHHG	RDSL E V L F Q G	240
241	PLAM D G F A G S	L D D S I S A A S T	S D V Q D R L S A L	E S R V Q Q Q E D E	I T V L K A A L A D	V L R R L A I S E D	300
301	H V A S V K K S V S	S K G Q P S P R A V	I P M S C I T N G S	G A N R K P S H T S	A V S I A G K E T L	S S A A K S G T E K	360
361	K K E K P Q G Q R E	K K E E S H S N D Q	S P Q I R A S P S P	Q P S S Q P L Q I H	R Q T P E S K N A T	P T K S I K R P S P	420
421	A E K S H N S W E N	S D D S R N K L S K	I P S T P K L I P K	V T K T A D K H K D	V I I N Q E G E Y I	K M F M R G R P I T	480
481	M F I P S D V D N Y	D D I R T E L P P E	K L K L E W A Y G Y	R G K D C R A N V Y	L L P T G E I V Y F	I A S V V V L F N Y	540
541	E E R T Q R H Y L G	H T D C V K C L A I	H P D K I R I A T G	Q I A G V D K D G R	P L Q P H V R V W D	S V T L S T L Q I I	600
601	G L G T F E R G V G	C L D F S K A D S G	V H L C V I D D S N	E H M L T V W D W Q	R K A K G A E I K T	T N E V V L A V E F	660
661	H P T D A N T I I T	C G K S H I F F W T	W S G N S L T R K Q	G I F G K Y E K P K	F V Q C L A F L G N	G D V L T G D S G G	720
721	V M L I W S K T T V	E P T P G K G P K V	Y R R K H Q E L Q A	Q M E M L Q S P E Y	K L S K L R T S T I	M T D Y N P N Y C F	780
781	A G K T S S I S D L	K E V P R K N I T L	I R G L G H G A F G	E V Y E G Q V S G M	P N D P S P L Q V A	V K T L P E V C S E	840
841	Q D E L D F L M E A	L I I S K L N H Q N	I V R C I G V S L Q	S L P R F I L L E L	M A G G D L K S F L	R E T R P R P S Q P	900
901	S S L A M L D L L H	V A R D I A C G C Q	Y L E E N H F I H R	D I A A R N C L L T	C P G P G R V A K I	G D F G M A R D I Y	960
961	R A S Y Y R K G G C	A M L P V K W M P P	E A F M E G I F T S	K T D T W S F G V L	L W E I F S L G Y M	P Y P S K S N Q E V	1020
1021	L E F V T S G G R M	D P P K N C P G P V	Y R I M T Q C W Q H	Q A P E D R P N F A I	I L E R I E Y C T Q	D P D V I N T A L P	1080
1081	I E Y G P L V E E E	E K V P V R P K D P	E G V P P L L V S Q	Q A K R E E E R S P	A A P P P L P T T S	S G K A A K K P T A	1140
1141	A E V S V R V P R G	P A V E G G H V N M	A F S Q S N P P S E	L H R V H G S R N K	P T S L W N P T Y G	S W F T E K P T K K	1200
1201	N N P I A K K E P H	E R G N L G L E G S	C T V P P N V A T G	R L P G A S L L L E	P S S L T A N M K E	V P L F R L R H F P	1260
1261	C G N V N Y G Y Q Q	Q G L P L E A A T A	P G A G H Y E D T I	L K S K N S M N Q P	G P		1320

1-218: GST Red: HIS6-tag Green: 3C blue: EML4ALK F: F1174L ALK mutation

EML4 ALK wt ¹ amino acid sequence							
1	MDGFAGSLDD	SISAASTSDV	QDRLSALESR	VQQQEDEITV	LKAALADVLR	RLAISEDHVA	60
61	SVKKSVS SKG	QPSRAVIPM	SCITNGSGAN	RKPSHTSAVS	IAGKETLSSA	AKSGTEKKKE	120
121	KPQGQREKKE	ESHSNDQSPQ	IRASPSQPQS	SQPLQIHRQT	PESKNATPTK	SIKRPSPAEK	180
181	SHNSWENSDD	SRNKLKIPS	TPKLIPKVTK	TADKHKDVII	NQEGEYIKMF	MRGRPITMFI	240
241	PSVDVNYDDI	RTELPPEK LK	LEWAYGYR GK	DCRANVYLLP	TGEIVYFIAS	VVLFNYEER	300
301	TQRHYLGHTD	CVKCLAIHPD	KIRIATGQIA	GVDKDGRPLQ	PHVRVWDSVT	LSTLQIIGLG	360
361	TFERGVGCLD	FSKADSGVHL	CVIDDSNEHM	LTVWDWQKKA	KGAEIKTTNE	VVLAVEFHPT	420
421	DANTIITCGK	SHIFFWTW SG	NSLTRKQGIF	GKYEKPKFVQ	CLAFLGNGDV	LTGDSGGVML	480
481	IWSKTTVEPT	PGKGPVYRR	KHQELQAMQM	ELQSPEYKLS	KLRTSTIMTD	YNPNYCFAGK	540
541	TSSISDLKEV	PRKNITLIRG	LGHGAFGEVY	EGQVSGMPND	PSPLQVAVKT	LPEVCSEQDE	600
601	LDFLMEALII	SKFNHQNIVR	CIGVSLQSLP	RFILLELMAG	GDLKSFLRET	RPRPSQPSSL	660
661	AMLDLLHVAR	DIACGCQYLE	ENHFIHRDIA	ARNCLLTCPG	PGRVAKIGDF	GMARDIYRAS	720
721	YYRKGGCAML	PVKWMPPEAF	MEGIFTSKTD	TWSFGVLLWE	IFSLGYMPYP	SKSNQEVLEF	780
781	VTSGGRMDPP	KNCPGPVYRI	MTQCWQHQPE	DRPNFAIILE	RIEYCTQDPD	VINTALPIEY	840
841	GPLVEEEEKV	PVRPKDPEGV	PPLLVSQQAK	REEERSPAAP	PPLPTTSSGK	AAKKPTAAEV	900
901	SVRVPRGPAV	EGGHVNMAFS	QSNPPSELHR	VHGSRNKPTS	LWNPTYGSWF	TEKPTKKNNP	960
961	IAKKEPHERG	NLGLG S C T V	PPNVATGRLP	GASL L L E P S S	LTANMKEVPL	FRLRHFPCGN	1020
1021	VNYGYQQOGL	PLEAATAPGA	GHYEDTILKS	KNSMNQPGP			1080

bold letters: expressed part of EML4 (blue) and ALK (green) **RED** letters: variant in Fusionprotein

¹NCBI/Protein accession number BAF73611.1