

ProQinase™ EPHB3

EPH receptor B3

Recombinant Human Active Protein Kinase

HGNC Symbol: EPHB3

Synonyms: Hek2, Tyro6, ETK2

Product No.: 0231-0000-1

Lot: 001

Description: Human Kinase, C-terminal fragment, amino acids Q₅₈₅-V₉₉₈ (as in [NCBI/Protein](#) entry NP_004434.1), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: EPHB3 Lot 001, has been verified by mass spectrometry LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 80148 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: This kinase was not activated by special procedures

Storage buffer: 50 mM TRIS-HCl pH 8.0, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20 % glycerol

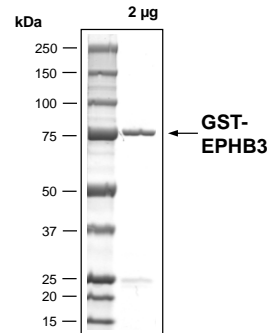
Storage temperature: -80°C

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!

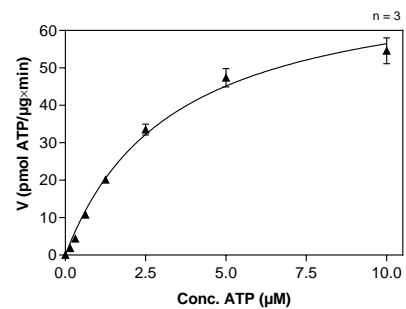
Protein concentration: 0.258 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:
Specific kinase activity (P_i transfer): 75 pmol/µg*min
ATP-K_M: 3.35 µM

EPHB3 Lot 001: Coomassie stain



EPHB3 Lot 001: Determination of V_{max} and K_M value for ATP



- 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: Poly(E/Y)_{4,1} 10 µg/ml
 - Kinase: 0.5 µg/ml

Assay technology:
Radiometric filter binding assay
MSFC membrane (96 well plate, Millipore)

Recombinant Proteins

Sequence information

GST-EPHB3 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEM	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPIQID	KYLKSSKYIA	WPLQGWAQATF	GGGDHPPKSD	PMGHHHHHGG	RRRASVAAGI	240
241	LVPRGSPGLD	GIYARGIQAS	MGARGRQCDG	YLQNSPLCMQ	RHGSDSEYTE	KLQYYIAPGM	300
301	KVYIDPFTYE	DPNEAVREFA	KEVDVSCVKI	EEVIGAGEFG	EVCGRRLKQP	GRREVFVAIK	360
361	TLKVGYTEERQ	RRDFLSEASI	MGQFDHPNII	RLEGVVTCSR	PVMILTEFME	NCALDSFLRL	420
421	NDGQFTVIQL	VGMLRGIAG	MKYLSEMYV	HRDLAARNIL	VNSNLVCKVS	DFGLSRFLED	480
481	DPSDPTYTSS	LGGKIPIRWT	APEAIAYRKF	TSASDVWSYG	IVMWEVMSYG	ERPYWDMNSQ	540
541	DVINAVEQDY	RLPPPMDCPT	ALHQLMLDCW	VRDRNLRPKF	SQIVNTLDKL	IRNAASLKVI	600
601	ASAQSGMSQP	LLDRTVPDYT	TFTTVGDWLD	AIKMGRYKES	FVSAGFASFD	LVAQMTAEDL	660
661	LRIGVTLAGH	QKILSSIQD	MRLQMNQTL	VQVKGEFQHT	GGRY		720

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site blue: EPHB3 fragment boxed: variation from RefSeq

EPHB3 wt ¹ Amino Acid Sequence									
1	MARARPPPPP	SPPPGLLPLL	PPLLLLPLLL	LPAGCRAL	EE TLM	DTKWVTS	ELAWTSHPES	60	
61	GWEEVSGYDE	AMNPIR	TYQV CNV	RESSQNN	WLRTGFI	WRR DVQR	VYVELK FT	VRDCNSIP	120
121	NIPGSCKETF	NLFYYE	ADSD VAS	ASSPFWM	ENPYVK	VDTI AP	DESFSRLD	AGRVNTK	VRS 180
181	FGPLSKAGFY	LAFQDQ	GACM SLIS	VRAFYK	KCASTT	AGFA LF	PETLTGAE	PTSLVI	APGT 240
241	CIPNAVEVSV	PLKLYC	NGDG EWM	VPVACT	CATGHE	PAAK ES	QCRPCPPG	SYKAKQ	GEGP 300
301	CLPCPPNSRT	TSPAAS	ICTC HNN	FYRADSD	SADSACT	TVP SPP	RGVISNV	NETSLI	LEWS 360
361	EPRDLGGRDD	LLYNVIC	KKC HG	AGGASACS	RCDDN	VEFVP RQ	LGLTERV	HISHLL	AHTR 420
421	YTFEVQAVNG	VSGKSPL	PPR YAA	VNITTNQ	AAPSEV	PTLR LH	SSSGSSLT	LSWAPP	ERP 480
481	GVILDYEMKY	FEKSEGI	AST VTS	QMNSVQL	DGLRP	DARYV VQ	VRARTVAG	YGQYSR	PAEF 540
541	ETTSEKSGA	QQLQEQL	PLI VGS	ATAGLVF	VVAVV	VIAIV CLR	KQRHGS	SEYTEKL	QY 600
601	IAPGMKVYID	PFTYED	PNEA VRE	FAKEIDV	SCVKIE	EVIG AGE	FGEVCRG	RLKQPG	RE 660
661	FVAIKTLKVG	YTERQRR	DFL SE	SIMQFD	HPNIIR	LEGV VT	KSRPVMIL	TEFMEN	CALD 720
721	SFLRLNDGQF	TVIQLV	GMLR GIA	AGMKYLS	EMNYV	HRDLA AR	NILVNSNL	VCKVSD	FGLS 780
781	RFLEDDPSDP	TYTSSL	GKPI PIR	WTAPEAI	AYRKFT	SASD VWS	YGIVMWE	VMSYGER	PYW 840
841	DMSNQDVINA	VEQDYRL	PPP MDC	PALHQL	MLDCW	VRDRN LR	PKFSQIVN	TLDKLI	RNAA 900
901	SLKVIASAQS	GMSQPL	LDRT VPD	YTTFTTV	GDWLD	AIKMG RY	KESFVSAG	FASFDL	VQA 960
961	TAEDLLRIGV	TLAGHQ	KKIL SSI	QDMRLQ	M NQ	TLFVQV			1020

blue: EPHB3 sequence expressed in recombinant protein Red: variant in recombinant protein

¹NCBI/Protein accession number NP_004434.1