

FAK aa2-1052

focal adhesion kinase

Recombinant Human Active Protein Kinase

HGNC Symbol: PTK2

Synonyms: FADK, FADK1, FAK1, FRNK, pp125FAK

Product No.: 0165-0000-3

Lot: 016

Description: Human FAK, amino acids A₂-H₁₀₅₂ (as in NCBI/Protein entry NP_722560.1), activated, N-terminal GST fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: FAK Lot 016, was confirmed as FAK by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 145,394 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: in-vitro activation by SRC

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: 0.249 µg/µl

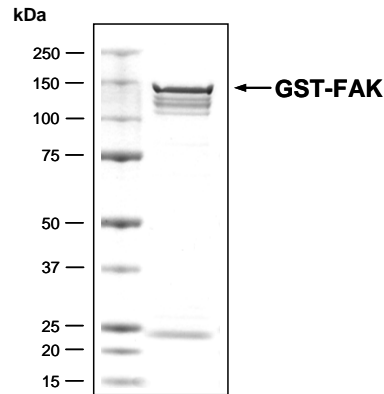
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

Specific kinase activity (P_i transfer): 14 pmol/µg × min

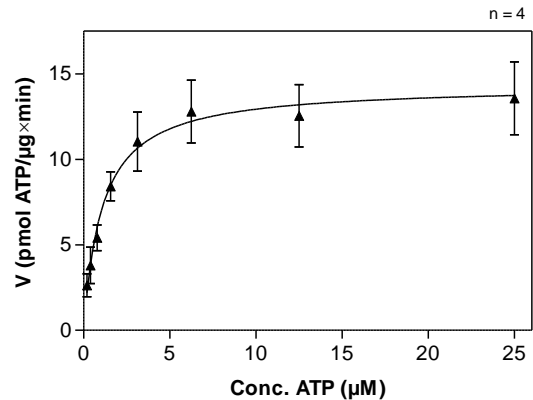
ATP-K_M: 1.1 µM

**FAK Lot 016:
Coomassie stain**



2.0 µg GST-FAK

**FAK Lot 016:
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: Poly(Glu/Tyr)_{4:1} 5 µg/ml
 - Kinase: 1 µg/ml
- Filter binding assay
 - MSFC membrane (Millipore)

Additional assay technology: FAK Lot 016

was also successfully tested by ProQinase 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



FAK

Product No.: 0165-0000-3

GST-FAK 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	LVPRGSAAAY	LDPNLNHTPN	240
241	SSTKTHLGTG	MERSPGAMER	VLKVFHYFES	NSEPTTWASI	IRHGATDVR	GIIQKIVDSH	300
301	KVKHVACYGF	RLSHLRSEEV	HWLHVDMGVS	SVREKYEALH	PPEEWKYELR	IRYLPKGFLN	360
361	QFTEDKPTLN	FFYQQVKSDY	MLEIADQVDQ	EIALKLGCLE	IRRSYWEMRG	NALEKKSNYE	420
421	VLEKDVGLKR	FFPKSLDSDV	KAKTLRKLIQ	QTFRQFANLN	REESILKFFE	ILSPVYRFDK	480
481	ECFKALGSS	WIISVELAIG	PEEGISYLT	KGCNPTHLD	FTQVQTIQYS	NSEDKDRKGM	540
541	LQLKIAGAPE	PLTVTAPSLT	IAENMADLID	GYCRLVNGTS	QSFIIRPQKE	GERALPSIPK	600
601	LANSEKQGMR	THAVSVSETD	DYAEIIDEED	TYTMPSTRDY	EIQRERIELG	RCIGEGQFGD	660
661	VHQGIYMSPE	NPALAVAIKT	CKNCTSDSVR	EKFLQEALTM	RQFDHPHIVK	LIGVITENPV	720
721	WIIMELCTLG	ELRSFLQVRK	YSLDLASLIL	YAYQLSTALA	YLESKRFVHR	DIAARNVLVS	780
781	SNDCVKLGDF	GLSRYMEDST	YKASKGKLP	IKWMAPESIN	FRFRTSASDV	WMFGVCMWEI	840
841	LMHGKVPFQG	VKNNDVIGRI	ENGERLPMPP	NCPPTLYSLM	TKCWAYDSPR	RPRFTELKAI	900
901	LSTILEEKA	QOEERMRMES	RRQATVSWDS	GGSDDEAPPK	SRPGYPSPRS	SEGFYPSQOH	960
961	MVQTNHYQVS	GYPGSHGITA	MAGSIYPGQA	SLLDQTDSDN	HRPQEIAMWQ	PNVEDSTVLD	1020
1021	LRGIGQVLP	HLMEERLIRQ	QOEMEEDQRW	LEKEERFLKP	DVRLSRGSID	REDGSLQGP	1080
1081	GNQHIYQVPG	KPDPAAPPK	PPRPGAPGHL	GSLASLSSPA	DSYNEGVLKQ	PQEISPPPTA	1140
1141	NLDRSNDKVY	ENVTGLVKAV	IEMSSKIQPA	PPEEYVPMVK	EVGLALRLL	ATVDETIPLL	1200
1201	PASTHREIEM	AQKLLNSDLG	ELINKMKLAQ	QYVMTSLQOE	YKQMLTAAH	ALAVDAKNLL	1260
1261	DVIDQARLKM	LGQTRPH					1320

1-218: GST **Pink**: Thrombin cleavage site **blue**: FAK **boxed**: variation from RefSeq

FAK wt ¹ Amino Acid Sequence							
1	MAAAYLDPNL	NHTPNSSTKT	HLGTGMERSP	GAMERVLKVF	HYFESNSEPT	TWASIIRHGD	60
61	ATDVRGIIQK	IVDSHKVKHV	ACYGFRLSHL	RSEEVHVLHV	DMGVSSVREK	YELAHPEEWE	120
121	KYELRIRYLP	KGFLNQFTED	KPTLNFFYQQ	VKSDYMLEIA	DQVDQEIALK	LGCLEIRRSY	180
181	WEMRGNALEK	KSNYEVLEKD	VGLKRFFPKS	LLDSVKAKTL	RKLIQQTFRQ	FANLNREESI	240
241	LKFFEILSPV	YRFDKECFKC	ALGSSWIIISV	ELAIGPEEGI	SYLTDKGCNP	THLADFTQVQ	300
301	TIQYSNSEDK	DRKGMQLKI	AGAPEPLTVT	APSLTIAENM	ADLIDGYCRL	VNGTSQSFI	360
361	RPQKEGERAL	PSIPKLANSE	KQGMRTHAVS	VSETDDYAEI	IDEEDTYTMP	STRDYEIQRE	420
421	RIELGRCIGE	GQFGDVHQGI	YMSPENPALA	VAIKTCKNCT	SDSVREKFLQ	EALTMRFQDH	480
481	PHIVKLIGVI	TENPVWIIME	LCTLGELRSF	LQVRKYSLDL	ASLILYAYQL	STALAYLESK	540
541	RFVHRDIAAR	NVLVSSNDCV	KLGDVGLSRY	MEDSTYYKAS	KGKLPKWMA	PESINFRRFT	600
601	SASDVWFMGV	CMWEILMHGV	KPFQGVKNND	VIGRIENGER	LPMPNPCPPT	LYSLMTKCWA	660
661	YDPSRRPRFT	ELKAQLSTIL	EEEKAQQEER	MRMESRRQAT	VSWDSGGSDE	APPKPSRPGY	720
721	PSPRSSEGFY	PSPQHMVQTN	HYQVSGYPGS	HGITAMAGSI	YPGQASLLDQ	TDSWNHRPQE	780
781	IAMWQPNVED	STVLDLRGIG	QVLPHTLMEE	RLIRQQQEME	EDQRWLEKEE	RFLKPDVRLS	840
841	RGSIDREDGS	LQGPIGNQHI	YQVPGKPDPA	APPKPPRPG	APGHLGSLAS	LSSPADSYNE	900
901	GVKLPQOEIS	PPPTANLDRS	NDKVYENVTG	LVKAVIEMSS	KIQPAPPEEY	VPMVKEVGLA	960
961	LRTLLATVDE	TIPLLPASTH	REIEMAQKLL	NSDLGELINK	MKLAQQYVMT	SLQOEYKQOM	1020
1021	LTAHALAVD	AKNLLDVIDQ	ARLKMVGQTR	PH			1080

blue: FAK sequence expressed in recombinant protein

¹NCBI/Protein accession number NP_722560.1