

## ProQinase™ VEGFR1

fms related receptor tyrosine kinase 1

Recombinant Human Active Protein Kinase

HGNC Symbol: FLT1

Synonyms: FLT, VEGF-R1

Product No.: 0097-0000-3

Lot: 011

**Description:** Human VEGFR1, C-terminal fragment, amino acids K<sub>784</sub>-I<sub>1338</sub> (as in [NCBI/Protein](#) entry NP\_002010.1), N-terminal GST-HIS<sub>6</sub> fusion protein with a Factor Xa cleavage site, expressed in Sf9 insect cells

**Product identity:** VEGFR1 LOT011, has been verified by mass spectrometry LC-ESI-MS/MS

**Theoretical MW**<sub>Fusion Protein</sub>: 89357 Da

**Expression host:** Sf9 insect cells

**Purification:** GST-Affinity Chromatography

**Activation:** in vitro auto activation

**Storage buffer:** 50 mM TRIS-HCl pH 8.0, 100 mM NaCl, 5 mM DTT, 4 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:** 0.16 µg/µl

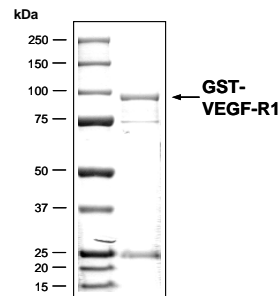
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Biochemical Parameters:**

Specific kinase activity (P<sub>i</sub> transfer): 151 pmol/µg\*min

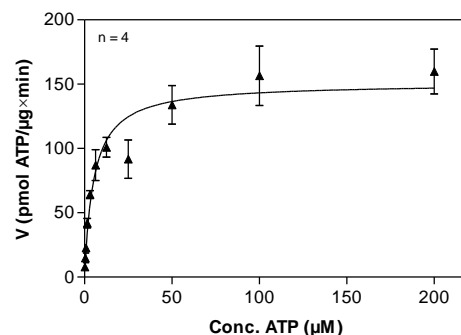
ATP-K<sub>M</sub>: 5.2 µM

### VEGFR1 LOT 011: Coomassie stain



2.0 µg GST-VEGF-R1

### VEGFR1 LOT 011: 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: Poly(Glu:Tyr)<sub>4:1</sub> 40 µg/ml
  - Kinase: 2 µg/ml

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

Sequence information

GST-VEGFR1 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	L <b>LE</b> GRGIL <b>KR</b>	<b>SSSEIKTDYL</b>	240
241	<b>SIIMDPDEVP</b>	<b>LDEQCERLPY</b>	<b>DASKWEFARE</b>	<b>RLKLGKSLGR</b>	<b>GAFGKVQAS</b>	<b>AFGIKKSPTC</b>	300
301	<b>RTVAVKMLKE</b>	<b>GATASEYKAL</b>	<b>MTELKILTHI</b>	<b>GHHLNVNLL</b>	<b>GACTKQGGPL</b>	<b>MVIVEYCKYG</b>	360
361	<b>NLSNYLKSQR</b>	<b>DLFFLNKDAE</b>	<b>LHMEPKKEKM</b>	<b>EPGLEQKPK</b>	<b>RLDVSTSES</b>	<b>FASSGFQEDK</b>	420
421	<b>SLSDVEEED</b>	<b>SDFGYKEPI</b>	<b>MEDLISYSFQ</b>	<b>VARGMEFLSS</b>	<b>RKCIHRDLAA</b>	<b>RNILLSENNV</b>	480
481	<b>VKICDFGLAR</b>	<b>DIYKNPDYVR</b>	<b>KGDTRLPLKW</b>	<b>MAPESIFDKI</b>	<b>YSTKSDVWSY</b>	<b>GVLLWEIFSL</b>	540
541	<b>GGSPYPGVQM</b>	<b>DEDFCSRLE</b>	<b>GMRMRAPEYS</b>	<b>TPEIQIMLD</b>	<b>CWHRDPKERP</b>	<b>RFAELVEKLG</b>	600
601	<b>DLLQANVQQD</b>	<b>GKDYIPINAI</b>	<b>LTGNSGFTYS</b>	<b>TPAFSEDFK</b>	<b>ESISAPKFN</b>	<b>GSSDDVRYVN</b>	660
661	<b>AFKFMSLERI</b>	<b>KTFEELLNA</b>	<b>TSMFDDYQD</b>	<b>SSTLLASPL</b>	<b>KRFTWTDSK</b>	<b>KASLKIDLRV</b>	720
721	<b>TSKSKESGLS</b>	<b>DVSRPSFCHS</b>	<b>SCGHVSEGKR</b>	<b>RFTYDHAELE</b>	<b>RKIACCSPPP</b>	<b>DYNSVLYST</b>	780
781	<b>PPI</b>						840

1-218: GST **Pink**: Factor Xa cleavage site **blue**: VEGFR1 fragment

VEGFR1 wt <sup>1</sup> Amino Acid Sequence							
1	MVSYWDTGVL	LCALLSCLLL	TGSSSGSKLK	DPESLKGTO	HIMQAGQTLH	LQCRGEAAHK	60
61	WSPPEMVSKE	SERLSITKSA	CGRNGKQFCS	TLTLNTAQAN	HTGFYSCYYL	AVPTSCKKET	120
121	ESAIYIFISD	TGRPFVEMYS	EIPEIIHMT	GRELVIKRV	TSPNITVTLK	KFPLDTLIPD	180
181	GKRIIWDNRK	GFIISNATYK	EIGLLTCEAT	VNGHLYKTY	LTHRQTNTII	DVQISTPRPV	240
241	KLLRGHTLV	NCTATTPNLT	RVQMTWSYD	EKNKRASVRR	RIDQNSHAN	IFYSVLTIDK	300
301	MQNKDKGLYT	CRVRSGPSFK	SVNTSVHIYD	KAFITVKHRK	QQVLETVAGK	RSYRLSMKVK	360
361	AFPSPEVWVL	KDGLPATEKS	ARYLTRGYSL	IIKDVTEEDA	GNYTILLSIK	QSNVFNKNTA	420
421	TLIVNVKQPI	YEKAVSSFPD	PALYPLGSRQ	ILTCTAYGIP	OPTIKWFVHP	CNNHSEARC	480
481	DFCSNNEESF	ILDADSNMGN	RIESITQRMA	IEGKNKMAS	TLVVADSRIS	GIYICIASNK	540
541	VGTVGRNIF	YITDVPNGFH	VNLEKMPTEG	EDLKLCTVN	KFLYRDVTWI	LLRVTNNRTM	600
601	HYSISKQKMA	ITKEHSITLN	LTIMNVSLQD	SGTYACRARN	VYTGEIILQK	KEITIRDQEA	660
661	PYLLRNLSDH	TVAISSSTTL	DCHANGVPEP	QITWFKNNHK	IQQEPGIILG	PGSSTLFIER	720
721	VTEEDEGVYH	CKATNQKGSV	ESSAYLTVQG	TSKSNLELI	TLTCTCVAAT	LFWLLLLLLI	780
781	RKMKRSSSEI	<b>KTDYLSIIMD</b>	<b>PDEVPLDEQC</b>	<b>ERLPYDASKW</b>	<b>EFARERLKL</b>	<b>KSLGRGAFGK</b>	840
841	<b>VVQASAFGIK</b>	<b>KSPTCRTVAV</b>	<b>KMLKEGATAS</b>	<b>EYKALMTELK</b>	<b>ILTHIGHHLN</b>	<b>VNLLGACTK</b>	900
901	<b>QGGPLMVIVE</b>	<b>YCKYGNLSNY</b>	<b>LKSKRDLFFL</b>	<b>NKDAALHMEP</b>	<b>KKEKMPEGLE</b>	<b>QGKPRLDSV</b>	960
961	<b>TSSESFASSG</b>	<b>FQEDKSLSDV</b>	<b>EEEESDGFY</b>	<b>KEPITMEDLI</b>	<b>SYSFQVARGM</b>	<b>EFLSSRKCIH</b>	1020
1021	<b>RDLAARNILL</b>	<b>SENNVVKICD</b>	<b>FGLARDIYKN</b>	<b>PDYVRKGDTR</b>	<b>LPLKWMAPES</b>	<b>IFDKIYSTKS</b>	1080
1081	<b>DVWSYGVLLW</b>	<b>EIFSLGGSPY</b>	<b>PGVQMEDDFC</b>	<b>SRLREGMRMR</b>	<b>APEYSTPEIY</b>	<b>QIMLDCWHRD</b>	1140
1141	<b>PKERPRFAEL</b>	<b>VEKLGDLLQA</b>	<b>NVQDGDYI</b>	<b>PINAILTGNS</b>	<b>GFTYSTPAFS</b>	<b>EDFFKESISA</b>	1200
1201	<b>PKFNSGSSDD</b>	<b>VRYVNAFKFM</b>	<b>SLEIKTFEE</b>	<b>LLPNATSMFD</b>	<b>DYQDSSTLL</b>	<b>ASPMLKRFTW</b>	1260
1261	<b>TDSKPKASLK</b>	<b>IDLRVTSKSK</b>	<b>ESGLSDVSRP</b>	<b>SFCHSSCGHV</b>	<b>SEGKRRFTYD</b>	<b>HAELERKIAC</b>	1320
1321	<b>CSPPPDYNSV</b>	<b>VLYSTPPI</b>					1380

**blue**: VEGFR1 sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](#) accession number NP\_002010.1