

ProQinase™ VEGFR2

vascular endothelial growth factor receptor 2

Recombinant Human Active Protein Kinase

HGNC Symbol: KDR

Synonyms: FLK1, VEGFR, CD309

Product No.: 0096-0000-1

Lot: 016

Description: Human VEGFR2, C-terminal fragment, amino acids D₈₀₇-V₁₃₅₆ (as in [NCBI/Protein](#) entry NP_002244.1), N-terminal GST fusion protein, expressed in Sf9 insect cells

Product identity: VEGFR2 Lot 016, was confirmed as VEGFR2 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 87,095 Da

Expression host: Sf9 insect 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

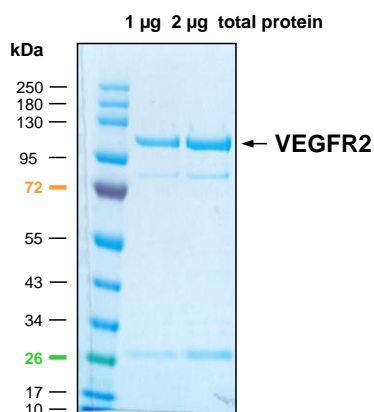
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.317 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

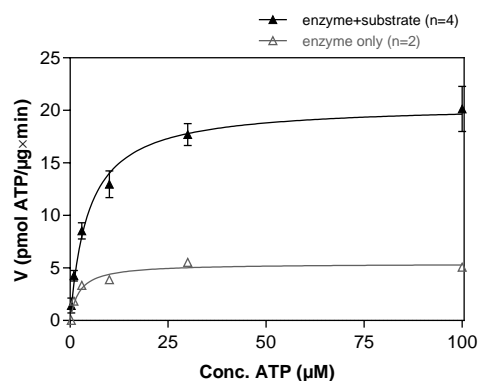
Biochemical Parameters:

Specific kinase activity (P_i transfer): 21 pmol/µg x min
ATP-K_M: 4.7 µM

VEGFR2 Lot 016: Coomassie stain



VEGFR2 Lot 016: 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(Glu:Tyr)_{4:1} 20 µg/ml
 - Kinase: 4 µg/ml
- Filter binding assay
MSFC membrane (Millipore)

Additional assay technology:

VEGFR2 Lot 016 was also successfully tested by Reaction Biology for the use with the ADP-Glo Kinase assay from Promega. ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

VEGFR2

Product No.: 0096-0000-1

Recombinant Proteins

GST-VEGF-R2 Recombinant Fusion Protein Amino Acid Sequence

1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGQATF	GGGDHPPKSD	PDELPLDEHC	ERLPYDASKW	240
241	EFPRDLRLKG	KPLGRGAFGQ	VIEADAFGID	KTATCRTVAV	KMLKEGATHS	EHRALMSELK	300
301	ILIHIGHHLN	VVNLGACTK	PGGPLMVIVE	FCKFGNLSTY	LRSKRNEFVP	YKTKGARFRQ	360
361	GKDYVGAIPV	DLKRRLDST	SSQSSASSGF	VEEKSLSDVE	EEEAPEDLYK	DFLTLEHLIC	420
421	YSFQVAKGME	FLASRCKIHR	DLAARNILLS	EKNVVKICDF	GLARDIYKDP	DYVRKGDARL	480
481	PLKWMAPETI	FDRVYTIQSD	VWSFGVLLWE	IFSLGASPYP	GVKIDEEFCR	RLKEGTRMRA	540
541	PDYTTPEMYQ	TMLDCWHGEP	SQRPTFSELV	EHLGNLLQAN	AQQDGDYIV	LPISSETLSME	600
601	EDSGLSLPTS	PVSCMEEEV	CDPKFHYDNT	AGISQYLQNS	KRKS RPVSVK	TFEDIPLLEP	660
661	EVKVIPDDNQ	TDSGMVLASE	ELKTLEDRTK	LSPSFGGMVP	SKSRESVASE	GSNQTSGYQS	720
721	GYHSDTDTT	VYSSEEAEEL	KLIEIGVQTG	STAQILQPDS	GTTLSSPPV		780

1-218: GST blue: VEGF-R2 fragment

VEGF-R2 wt¹ Amino Acid Sequence

1	MQSKVLLAVA	LWLCVETRAA	SVGLPSVSLD	LPRLSIQKDI	LTIKANTTLQ	ITCRGQRDL	60
61	WLWPNNQSGS	EQRVEVTECS	DGLFCKTLTI	PKVIGNDTGA	YKCFYRETDL	ASVIYVYVQD	120
121	YRSPFIASVS	DQHGCVVYITE	NKNKTVVIPC	LGSISNLNVS	LCARYPEKRF	VPDGNRISWD	180
181	SKKGFTIPSY	MISYAGMVFC	EAKINDESYQ	SIMYIVVVVG	YRIYDVVLSF	SHGIELSVGE	240
241	KLVLNCTART	ELNVGIDFNW	EYPSSKHQHK	KLVNRLDKTQ	SGSEMKKFLS	TLTIDGVTRS	300
301	DQGLYTCAAS	SGLMTKKNST	FVRVHEKPFV	AFGSGMESLV	EATVGERVRI	PAKYLGYPPP	360
361	EIKWYKNGIP	LESNHTIKAG	HVLTIMEVSE	RDTGNYTVIL	TNPISKEKQS	HVVSLVVYVP	420
421	PQIGEKSLIS	PVDSYQYGT	QTLTCTVYAI	PPPHIHWHY	QLEEECANEP	SQAVSVTNPY	480
481	PCEEWRVED	FQGGNKIEVN	KNQFALIEGK	NKTVSTLVIQ	AANVSALYKC	EAVNKVGRGE	540
541	RVISFHVTRG	PEITLQPD	PTEQESVSLW	CTADRSTFEN	LTWYKLGQP	LPIHVGELPT	600
601	PVCKNLDTLW	KLNATMFSNS	TNDILIMELK	NASLQDQGDY	VCLAQDRRTK	KRHCVVRQLT	660
661	VLERVAPTIT	GNLENQTTSI	GESIEVSC	SGNPPPQIMW	FKDNETLVED	SGIVLKDGNR	720
721	NLTIRVRKE	DEGLYTCQAC	SVLGC	AKVEA	FFIIEGAQEK	TNLEIIILVG	780
781	LLVIIILRTVK	RANGGELKTG	YLSIVMDPDE	LPLDEHCERL	PYDASKWEFP	RDRLKLGKPL	840
841	GRGAFGQVIE	ADAFGIDKTA	TCRTVAVKML	KEGATHSEHR	ALMSELKILI	HIGHHLNVN	900
901	LLGACTKPGG	PLMVIVEFCK	FGNLSTYLRS	KRNEFVPYKT	KGARFRQGD	YVGAIPVDLK	960
961	RRLDSITSSQ	SSASSGFVEE	KSLSDVEEEE	APEDLYKDFL	TLEHLICYSF	QVAKGMFLA	1020
1021	SRKCIHRDLA	ARNILLSEKN	VVKICDFGLA	RDIYKDPDYV	RKGDARLPLK	WMAPETIFDR	1080
1081	VYTIQSDVWS	FGVLLWEIFS	LGASPYPGVK	IDEEFCRRLK	EGTRMRAPDY	TPPEMYQTML	1140
1141	DDTHGEP	PTFSELVEHL	GNLLQANAQQ	DGKDYIVLPI	SETLSMEEDS	GLSLPTSPVS	1200
1201	CMEEEEVCDP	KFHYDNTAGI	SQYLQNSKRK	SRPVSVKTFE	DIPLEEPEVK	VIPDDNQ	1260
1261	GMVLASEELK	TLEDRTKLSP	SFGGMVPSKS	RESVASEGSN	QTSYQSGYH	SDDTDTTVYS	1320
1321	SEEAEELKLI	EIGVQTGSTA	QILQPD	SGTT	LSSPPV		1380

blue: VEGF-R2 sequence expressed in recombinant protein

¹NCBI/Protein accession number NP_002244.1