

ProQinase™ CCDC6-RET

CCDC6-RET fusion protein variant a

Recombinant Human Active Protein Kinase

HGNC Symbol: n/a

Synonyms: CCDC6-RETA

Product No.: 1417-0000-1

Lot: 003

Description: Human CCDC6-RET fusion protein, amino acids Met₁-Ser₅₀₃ (as in [NCBI/Protein](#) entry BAM36435.1), N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: CCDC6-RET Lot 003, was confirmed as a CCDC6-RET by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 84,394 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: 0.178 µg/µl

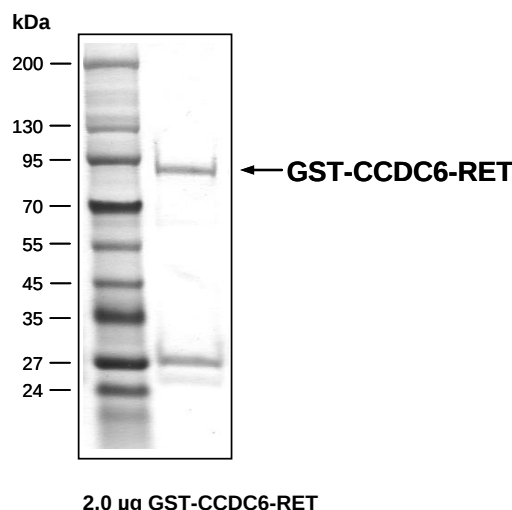
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

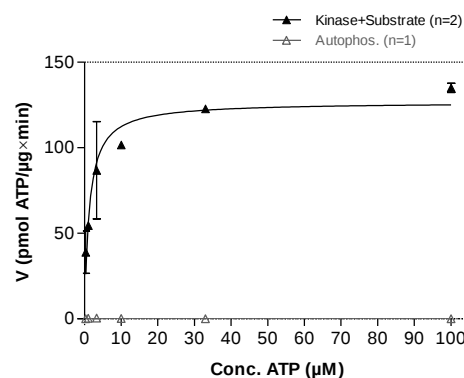
Specific kinase activity (P_i transfer): 127 pmol/µg × min

ATP-K_M: 1.3 µM

CCDC6-RET Lot 003: Coomassie stain



CCDC6-RET Lot 003: 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: TRK-C derived peptide, 40 µg/ml
Kinase: 1.0 µg/ml
- Filter binding assay
MSPH membrane (Millipore)

Additional assay technology:

CCDC6-RET Lot 003 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

ProQinase™ CCDC6-RET

Product No.: 1417-0000-1

GST-CCDC6-RET Recombinant Fusion Protein Amino Acid Sequence

1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLIERDEGDK	WRNKKFELGL	EFNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPEML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVLFGG	240
241	PLAMGARGRM	ADSASESDTD	GAGGNSSSSA	AMQSSCSSTS	GGGGGGGGGG	GGGKSGGIVI	300
301	SPFRLEELTN	RLASLQENK	VLKIELETYK	LKCKALQEEN	RDLRKASVTI	EDPKWEFPRK	360
361	NLVLGKTLGE	GEFGKVVKAT	AFHLKGRAGY	TTVAVKMLKE	NASPSELRLD	LSEFNVLKQV	420
421	NHPHVIKLYG	ACSQDGPLLL	IVEYAKYGS	RGFLRESRKV	GPGYLGSGGS	RNSSSLDHPD	480
481	ERALTMGDLI	SFAWQISQGM	QYLAEMKLVH	RDLAARNILV	AEGRKMKISD	FGLSRDVYEE	540
541	DSYVKRSQGR	IPVKWMAIES	LFDHIYTTQS	DVWSFGVLLW	EIVTLGGNPY	PGIPPERLFN	600
601	LLKTGHRMER	PDNCSEEMYR	LMLQCWKQEP	DKRPVFADIS	KDLEKMMVKR	RDYDLAAS	660
661	PSDSLIIYDDG	LSEETPLVD	CNNAPLPRAL	PSTWIENKLY	GMSDPNWPGE	SPVPLTRADG	720
721	TNTGFPRYPN	DSVYANWMLS	PSAAKLMDTF	DS			780

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: CCDC6-RET

CCDC6-RET wt¹ Amino Acid Sequence

1	MADSASESDT	DGAGGNSSSS	AAMQSSCSST	SGGGGGGGGG	GGGKSGGIV	ISPFRLLELT	60
61	NRLASLQEN	KVLKIELETY	KLKCKALQEE	NRDLRKASVT	IEDPKWEFPR	KNLVLGKTLG	120
121	EGEFGKVVK	TAFHLKGRAG	YTTVAVKMLK	ENASPSELRLD	LLSEFNVLKQ	VNHPHVIKLY	180
181	GACSDGPLL	LIVEYAKYGS	LRGFLRESRK	VGPGYLGSGG	SRNSSSLDHP	DERALTMGDL	240
241	ISFAWQISQG	MQYLAEMKLV	HRDLAARNIL	VAEGRKMKIS	DFGLSRDVYE	EDSYVKRSQG	300
301	RIPVKWMAIE	SLFDHIYTTQ	SDVWSFGVLL	WEIVTLGGNP	YPGIPPERLF	NLLKTGHRME	360
361	RPDNCSEEMY	RLMLQCWKQE	PDKRPVFADI	SKDLEKMMVK	RRDYDLAAS	TPSDSLIYDD	420
421	GLSEETPLV	DCNNAPLPRA	LPSTWIENKL	YGMSDPNWPGE	ESPVPLTRAD	GTNTGFPRYP	480
481	NDSVYANWML	SPSAAKLMDT	FDS				540

blue: CCDC6-RET sequence expressed in recombinant protein E₁₀₂: End of CCDC6 / start of RET ORF

¹NCBI/Protein accession number BAM36435.1