

FGF-R3 G697C

fibroblast growth factor receptor 3

Recombinant Human Active Protein Kinase

HGNC Symbol: FGFR3

Synonyms: ACH, CD333, CEK2, FGFR-3, HSVGFR3EX, JTK4

Product No.: 1071-0000-1

Lot: 001

Description: Human FGF-R3, C-terminal fragment, amino acids R₃₉₇-T₈₀₆ (as in NCBI/Protein entry NP_000133.1) with a G697C mutation, N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: FGF-R3 G697C, Lot 001, was confirmed as FGF-R3 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 73,980 Da

Expression: Baculovirus infected Sf9 cells

Purification: GST-Affinity Chromatography

Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20% glycerol

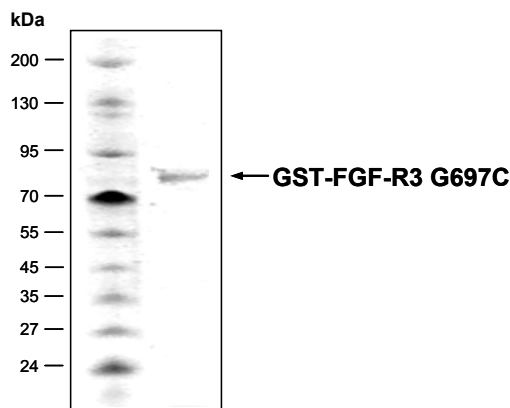
Storage temperature: -80°C
Avoid repeated freeze-thaw cycles!

Protein concentration: 0.225 µg/µl (Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

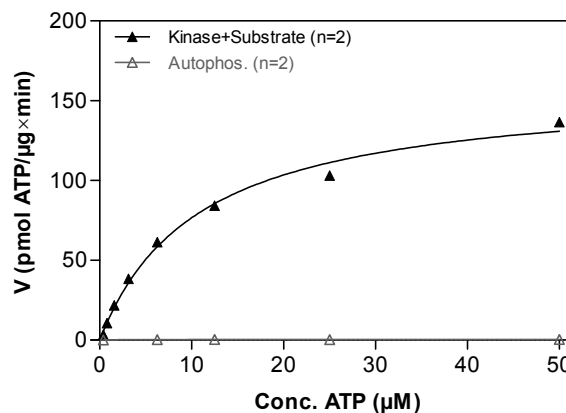
Specific activity: 159 pmol/µg×min
ATP-K_M: 11 µM

**FGF-R3 G697C Lot 001:
Coomassie stain**



2.0 µg GST-FGF-R3 G697C


**FGF-R3 G697C Lot 001:
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: TRK-C-derived Peptide (R11-VYSTDYRFLNPS), 40 µg / ml
 - FGF-R3 G697C: 1.0 µg / ml
- Filter binding assay
 - MSPH membrane (Millipore)

Additional assay technology: FGF-R3 G697C Lot 001

was also successfully tested by ProQinase for the use with the ADP-Glo™ Kinase assay from 

This product is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.

FGF-R3 G697C

Product No.: 1071-0000-1

FGF-R3 G697C Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQ SMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHHG	RDSLEVL FQG	240
241	PLAMVRLRSP	PKKGLGSPTV	HKISRFP LKR	QVSLESNASM	SSNTPLVRIA	RLSSGEGPTL	300
301	ANVSELELPA	DPKWELSRAR	LTLGKPLGEG	CFGQVVM AEA	IGIDKDRAAK	PVTVAVKMLK	360
361	DDATDKDLS	LVSEMEMMKM	IGKHKNIINL	LGACTQGGPL	YVLVEYAAKG	NLREFLRARR	420
421	PPGLDYSFDT	CKPPEEQ LTF	KDLVSCAYQV	ARGMEYLA SQ	KCIHRDLAAR	NVLVTEDNVM	480
481	KIADFGLAR	VHNL DYYKKT	TNGRLPVKWM	APEALFDRVY	THQSDVWSFG	VLLWEIFTLG	540
541	GSPYPCIPVE	ELFKLLKEGH	RMDK PANCTH	DLYMIMRECW	HAAPSQRPTF	KQLVEDLDRV	600
600	LTVTSTDEYL	DLSAPFEQYS	PGGQDTPSS	SSGDDSVFAH	DLLPPAPSS	GGSR T	660

1-218: GST **Red:** HIS6-tag **Green:** 3C **blue:** FGF-R3 fragment **C:** G697C mutation

FGF-R3 wt ¹ amino acid sequence							
1	MGAPACALAL	CVAVAIVAGA	SSESLGTEQR	VVGRAAEVPG	PEPGQQEQLV	FGSGDAVELS	60
61	CPPPGGGPMG	PTVWVKDGTG	LVPSE RVLVG	PQRLQVLNAS	HEDSGAYSCR	QRLTQRVLCH	120
121	FSVRVTDAPS	SGDDEDGEDE	AEDTGVD TGA	PYWTRPERMD	KLLAVPAAN	TVRFRCPAAG	180
181	NPTPSISWLK	NGREFRGEHR	IGGIKLRHQQ	WSLVMESVVP	SDRGNYTCVV	ENKFGSIRQT	240
241	YTLDV LERSP	HRPILQAGLP	ANQTAVLGSD	VEFHCKVYSD	AQPHIQWLKH	VEVNGSKVGP	300
301	DGTPYVTVLK	TAGANTTDKE	LEVLSLHNVT	FEDAGEYTCL	AGNSIGFSHH	SAWLVVLP AE	360
361	EELVEADEAG	SVYAGILSYG	VGFFLFILVV	AAVTLCRLRS	PPKKGLGSPT	VHKISRFP LK	420
421	RQVSLESNAS	MSSNTPLVRI	ARLSSGEGPT	LANVSELELP	ADPKWELSR A	RLTLGKPLGE	480
481	GCFGQVMAE	AIGIDKDRAA	KPVTVAVKML	KDDATDKDLS	DLVSEMEMMK	MIGKHKNIIN	540
541	LLGACTQGGP	LYVLVEYAAK	GNLREFLRAR	RPPGLDYSFD	TCKPPEEQ LTF	FKDLVSCAYQ	600
600	VARGMEYLA S	QKCIHRDLAA	RNVLVTEDNV	MKIADFGLAR	DVHNL DYYKKT	TTNGRLPVKW	660
661	MAPEALFDRV	YTHQSDVWSF	GVLLEIFTL	GGSPYCGIPV	EELFKLLKEG	HRMDK PANCT	720
721	HDLYMIMREC	WHAAPSQRPT	FKQLVEDLDR	VLTVTSTDEY	LDLSAPFEQY	SPGGQDTPSS	780
781	SSGDDSVFAH	HDLLPPAPPS	SGGSRT				840

blue: fragment of FGF-R3 expressed in fusionprotein **Red:** variant in fusionprotein

¹NCBI/Protein accession number NP_000133.1