

## FGF-R3 G697C

fibroblast growth factor receptor 3

Recombinant Human Active Protein Kinase

**HGNC Symbol:** FGFR3

**Synonyms:** ACH, CD333, CEK2, FGFR-3, HSFGR3EX, JTK4

**Product No.:** 1071-0000-1

**Lot:** 001

**Description:** Human FGF-R3, C-terminal fragment, amino acids R<sub>397</sub>-T<sub>806</sub> (as in [NCBI/Protein](#) entry NP\_000133.1) with a G697C mutation, N-terminal GST-HIS<sub>6</sub> 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<sub>Fusion Protein</sub>:** 73,980 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:** 0.225 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

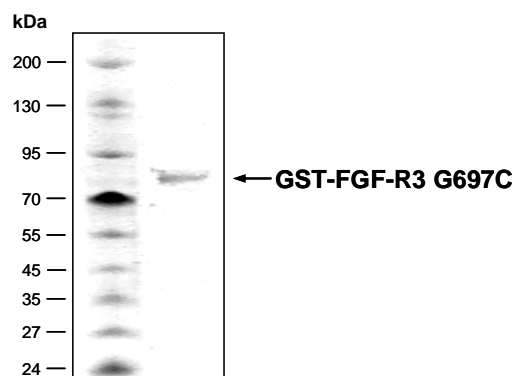
### Biochemical Parameters:

Specific kinase activity (P<sub>i</sub> transfer): 159 pmol/µg × min  
ATP-K<sub>M</sub>: 11 µM

### Additional assay technology:

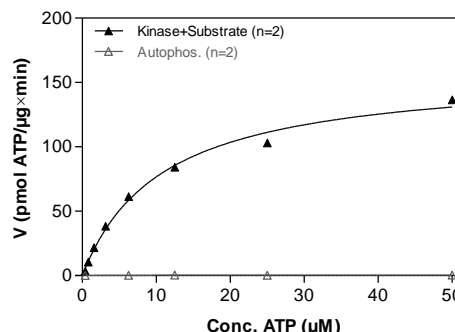
FGF-R3 G697C Lot 001 was also successfully tested by ProQinase for the use with the ADP-Glo™ Kinase assay from Promega  
ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

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



2.0 µg GST-FGF-R3 G697C

### FGF-R3 G697C Lot 001: 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: TRK-C derived peptide 40 µg/ml
  - Kinase: 1 µg/ml
- Filter binding assay  
MSPH membrane (Millipore)

## FGF-R3 G697C

Product No.: 1071-0000-1

GST-FGF-R3 G697C 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	PMG <b>HHHHHG</b>	RDS <b>LEVLFQG</b>	240
241	<b>PLAMVRLRSP</b>	<b>PKKGLGSPTV</b>	<b>HKISRFP LKR</b>	<b>QVSLESNASM</b>	<b>SSNTPLVRIA</b>	<b>RLSSGEGPTL</b>	300
301	<b>ANVSELELPA</b>	<b>DPKWELSRAR</b>	<b>LTLGKPLGEG</b>	<b>CFGQVMAEA</b>	<b>IGIDKDRAAK</b>	<b>PVTVAVKMLK</b>	360
361	<b>DDATDKDLS</b>	<b>LVSEMEMMKM</b>	<b>IGKHNIINL</b>	<b>LGACTQGGPL</b>	<b>YVLVEYAAKG</b>	<b>NLREFLRARR</b>	420
421	<b>PPGLDYSFDT</b>	<b>CKPPEEQLTF</b>	<b>KDLVSCAYQV</b>	<b>ARGMEYLASQ</b>	<b>KCIHRDLAAR</b>	<b>NVLVTEDNVM</b>	480
481	<b>KIADFG LARD</b>	<b>VHNLDYYKKT</b>	<b>TNGRLPVKWM</b>	<b>APEALFDRVY</b>	<b>THQSDVWSFG</b>	<b>VLLWEIFTLG</b>	540
541	<b>GSPYPCIPVE</b>	<b>ELFKLLKEGH</b>	<b>RMDKPANCTH</b>	<b>DLYMIMRECW</b>	<b>HAAPSQRPTF</b>	<b>KQLVEDLDRV</b>	600
600	<b>LTVTSTDEYL</b>	<b>DLSAPFEQYS</b>	<b>PGGQDTPSSS</b>	<b>SSGDDSVFAH</b>	<b>DLLPPAPPS</b>	<b>GGSR T</b>	660

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

FGF-R3 wt <sup>1</sup> Amino Acid Sequence							
1	MGAPACALAL	CVAVAIVAGA	SSESLGTEQR	VVGRAAEVPG	PEPGQEQLV	FGSGDAVELS	60
61	CPPPGGGPMG	PTVWVKDGTG	LVPSEVLVVG	PQRLQVLNAS	HEDSGAYSCR	QRLTQRVLCH	120
121	FSVRVTDAPS	SGDDEDGEDE	AEDTGVDTGA	PYWTRPERMD	KKLLAVPAAN	TVRFRCPAAG	180
181	NPTPSISWLK	NGREFRGEHR	IGGIKLRHQQ	WSLVMESVVP	SDRGNYTCVV	ENKFGSIRQT	240
241	YTLVDLERSP	HRPILQAGLP	ANQTAVLGSD	VEFHCKVYSD	AQPHIQWLKH	VEVNGSKVGP	300
301	DGTPYVTVLK	TAGANTTDEK	LEVLSLHNVT	FEDAGEYTCL	AGNSIGFSHH	SAWLVLPAE	360
361	EELVEADEAG	SVYAGILSYG	VGFFLFILVV	AAVTLCRLRS	<b>PPKKGLGSPT</b>	<b>VHKISRFP LK</b>	420
421	<b>RQVSLESNAS</b>	<b>MSSNTPLVRI</b>	<b>ARLSSGEGPT</b>	<b>LANVSELELP</b>	<b>ADPKWELSRA</b>	<b>RLTLGKPLGE</b>	480
481	<b>GCFGQVMAE</b>	<b>AIGIDKDRAA</b>	<b>KPVTVAVKML</b>	<b>KDDATDKDLS</b>	<b>DLVSEMEMMK</b>	<b>MIGKHNIIN</b>	540
541	<b>LLGACTQGGP</b>	<b>LYVLVEYAAK</b>	<b>GNLREFLRAR</b>	<b>RPPGLDYSFD</b>	<b>TCKPPEEQLT</b>	<b>FKDLVSCAYQ</b>	600
600	<b>VARGMEYLAS</b>	<b>QKCIHRDLAA</b>	<b>RNVLVTEDNV</b>	<b>MKIADFG LAR</b>	<b>DVHNLDYYK</b>	<b>TTNGRLPVKW</b>	660
661	<b>MAPEALFDRV</b>	<b>YTHQSDVWSF</b>	<b>GVLLWEIFTL</b>	<b>GGSPYPCIPV</b>	<b>EELFKLLKEG</b>	<b>HRMDKPANCT</b>	720
721	<b>HDLYMIMREC</b>	<b>WHAAPSQRPT</b>	<b>FKQLVEDLDR</b>	<b>VTVTSTDEY</b>	<b>LDLSAPFEQY</b>	<b>SPGGQDTPSS</b>	780
781	<b>SSSGDDSVFA</b>	<b>HDLLPPAPPS</b>	<b>SGGSRT</b>				840

**blue**: FGF-R3 sequence expressed in recombinant protein **Red**: variant in recombinant protein

<sup>1</sup>[NCBI/Protein](#) accession number NP\_000133.1

This product was manufactured at ProQinase in Freiburg, Germany, and 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.