

## ProQinase™ FGF-R3 K650M

fibroblast growth factor receptor 3

Recombinant Human Active Protein Kinase

HGNC Symbol: FGFR3

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

**Product No.:** 1069-0000-1

**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 K650M mutation, N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

**Product identity:** FGF-R3 K650M, Lot 001, was confirmed as FGF-R3 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>Fusion Protein</sub>:** 73,937 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

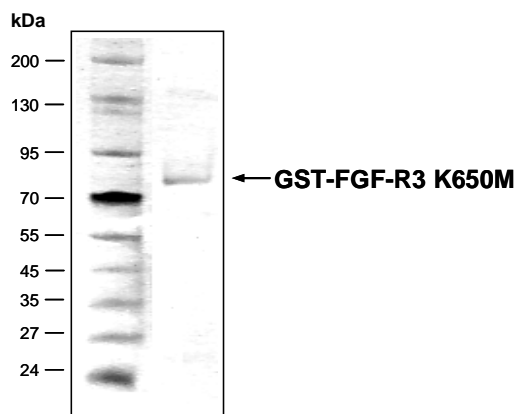
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

**Protein concentration:** 0.186 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Biochemical Parameters:**

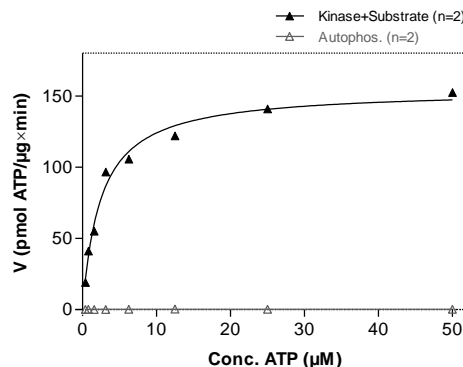
Specific kinase activity (P<sub>i</sub> transfer): 154 pmol/µg × min  
ATP-K<sub>M</sub>: 2.5 µM

**FGF-R3 K650M Lot 001:  
Coomassie stain**



2.0 µg GST-FGF-R3 K650M

**FGF-R3 K650M Lot 001:  
Determination of V<sub>max</sub> and K<sub>M</sub> value for ATP**



**Determination of K<sub>M</sub> value & Specific activity:**

- 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)

**Additional assay technology:**

FGF-R3 K650M Lot 001 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

This product was manufactured at Reaction Biology in Freiburg, Germany, and is for in vitro research use only, not for use in humans or animals.  
© European Union, 2020. Material may not be reproduced or distributed without written permission from Reaction Biology Europe GmbH

## ProQinase™ FGF-R3 K650M

Product No.: 01069-0000-1

| GST-FGF-R3 K650M Recombinant Fusion Protein Amino Acid Sequence |                   |                   |                    |                   |                   |                    |     |
|---|-------------------|-------------------|--------------------|-------------------|-------------------|--------------------|-----|
| 1   | MSPILGYWKI        | KGLVQPTRL         | LEYLEEKYEE         | HLYERDEGDK        | WRNKKFELGL        | EFPNLPYYID         | 60  |
| 61  | GDVKLTQSMA        | IIRYIADKHN        | MLGGCPKERA         | EISMLEGAVL        | DIRYGVSRIA        | YSKDFETLKV         | 120 |
| 121   | DFLSKLPPEML       | KMFEDRLCHK        | TYLNGDHVTH         | PDFMLYDALD        | VVLYMDPMCL        | DAFPKLVCFK         | 180 |
| 181   | KRIEAIPOID        | 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>LTGKPLGEG</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>YVLVEYAAK</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>KIADFGLARD</b> | <b>VHNLDYYKMT</b> | <b>TNGRLPVKWM</b>  | <b>APEALFDRVY</b> | <b>THQSDVWSFG</b> | <b>VLLWEIFTLG</b>  | 540 |
| 541   | <b>GSPYPGIPVE</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>DLLPPAPSSS</b> | <b>GGSR</b>        | 660 |

1-218: GST **Red**: HIS6-tag **Green**: 3C cleavage site **blue**: FGF-R3 fragment **boxed**: K650M

| 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>MKIADFGLAR</b> | <b>DVHNLDYYK</b>  | <b>TTNGRLPVKW</b>  | 660 |
| 661  | <b>MAPEALFDRV</b> | <b>YTHQSDVWSF</b> | <b>GVLLWEIFTL</b> | <b>GGSPYPGIPV</b> | <b>EELFKLLKEG</b> | <b>HRMDKPANCT</b>  | 720 |
| 721  | <b>HDLYMIMREC</b> | <b>WHAAPSQRPT</b> | <b>FKQLVEDLDR</b> | <b>VLTVTSTDEY</b> | <b>LDLSAPFEQY</b> | <b>SPGGQDTPSS</b>  | 780 |
| 781  | <b>SSSGDDSVFA</b> | <b>HDLLPPAPSS</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