

FGF-R3 wt

fibroblast growth factor receptor 3

Recombinant Human Active Protein Kinase

HGNC Symbol: FGFR3

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

Product No.: 1068-0000-1

Lot: 005

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

Product identity: FGF-R3 wt, Lot 005, was confirmed as FGF-R3 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 73,934 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: 0.302 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

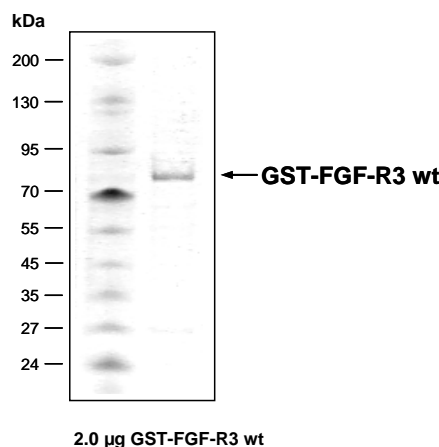
Biochemical Parameters:

Specific kinase activity (P_i transfer): 199 pmol/µg × min
ATP-K_M: 6.1 µM

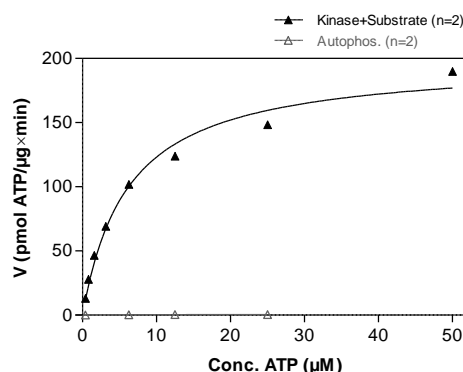
Additional assay technology:

FGF-R3 wt Lot 005 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 wt Lot 005: Coomassie stain



FGF-R3 wt Lot 005: 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, 40 µg/ml
 - Kinase: 1 µg/ml
- Filter binding assay
MSPH membrane (Millipore)

FGF-R3 wt

Product No.: 1068-0000-1

GST-FGF-R3 wt Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMG HHHHHG	RDS LEVLFCG	240
241	PLAMVRLRSP	PKKGLGSPTV	HKISRFP LKR	QVSLESNASM	SSNTPLVRIA	RLSSGEGPTL	300
301	ANVSELELPA	DPKWELSRAR	LTLGKPLGEG	CFGQVMAEA	IGIDKDRAAK	PVTVAVKMLK	360
361	DDATDKDLS	LVSEMEMMKM	IGKHNIINL	LGACTQGGPL	YVLVEYAAKG	NLREFLRARR	420
421	PPGLDYSFDT	CKPPEEQLTF	KDLVSCAYQV	ARGMEYLASQ	KCIHRDLAAR	NVLVTEDNVM	480
481	KIADFGLARD	VHNL DYYKKT	TNGRLPVKWM	APEALFDRVY	THQSDVWSFG	VLLWEIFTLG	540
541	GSPYPGIPVE	ELFKLLKEGH	RMDKPANCTH	DLYMIMRECW	HAAPSQRPTF	KQLVEDLDRV	600
600	LTVTSTDEYL	DLSAPFEQYS	PGGQDTPSSS	SSGDDSVFAH	DLLPPAPPS	GGSR T	660

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

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

blue: FGF-R3 sequence expressed in recombinant protein

¹[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.