

## FER

FER tyrosine kinase

Recombinant Human Active Protein Kinase

HGNC Symbol: FER

Synonyms: TYK3, c-FER, p94-FER, PPP1R74

Product No.: 0533-0000-1

Lot: 002

**Description:** Human FER, C-terminal fragment, amino acids K<sub>541</sub>-T<sub>822</sub> (as in [NCBI/Protein](#) entry NP\_005237.1), N-terminal GST-HIS<sub>6</sub> fusion protein with a Thrombin cleavage site, expressed in insect cells

**Product identity:** FER, Lot 002, was confirmed as FER by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW**<sub>Fusion Protein</sub>: 61,733 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, 4 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:** 0.183 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

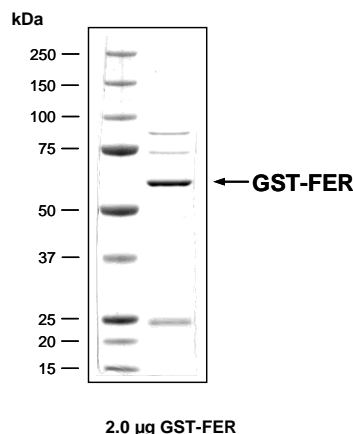
### Biochemical Parameters:

Specific kinase activity (P<sub>i</sub> transfer): 298 pmol/µg × min  
ATP-K<sub>M</sub>: 4.7 µM

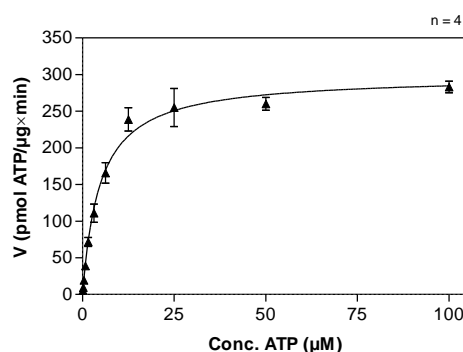
### Additional assay technology:

FER Lot 002 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

### FER Lot 002: Coomassie stain



### FER Lot 002: 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: Poly(Glu, Tyr)<sub>4:1</sub> 10 µg/ml  
Kinase: 4 µg/ml
- Filter binding assay  
MSFC membrane (Millipore)

# FER

Product No.: 0533-0000-1

| GST-FER 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  | KRIEAIPIQID | KYLKSSKYIA | WPLQGWQATF | GGGDHPPKSD | PMGHHHHHG  | RRRASVAAGI  | 240 |
| 241  | LVPRGSPGLD  | GICSRNSKSG | VVLLNPIPKD | KKWILSHEDV | ILGELLGKGN | FGEVYKGTLK  | 300 |
| 301  | DKTSVAVKTC  | KEDLPQELKI | KFLQEAAILK | QYDHPNIVKL | IGVCTQRQPV | YIIMELVSGG  | 360 |
| 361  | DFLTFLRRRK  | DELKLNKQV  | FSLDAAAGML | YLESKNCIHR | DLAARNCLVG | ENNVLKISDF  | 420 |
| 421  | GMSRQEDGGV  | YSSSGLKQIP | IKWTAPEALN | YGRYSSESDV | WSFGILLWET | FSLGVCPPYPG | 480 |
| 481  | MTNQAREQV   | ERGYRMSAPQ | HCPEDISKIM | MKCWDYKPEN | RPKFSELQKE | LTIIKRKLT   | 540 |

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site blue: FER fragment

| FER wt <sup>1</sup> Amino Acid Sequence |            |            |            |            |             |            |     |
|---|------------|------------|------------|------------|-------------|------------|-----|
| 1                                       | MGFGSDLKNS | HEAVLKLQDW | ELRLLETVKK | FMALRIKSDK | EYASTLQNL   | NQVDKESTVQ | 60  |
| 61                                      | MNYVSNVSKS | WLLMIQQTEQ | LSRIMKTHAE | DLNSGPLHRL | TMMIKDKQQV  | KKSYIGVHQQ | 120 |
| 121                                     | IEAEMIKVTK | TELEKLKCSY | RQLIKEMNSA | KEYKKEALAK | GKETEKAKER  | YDKATMKLHM | 180 |
| 181                                     | LHNQYVLALK | GAQLHQNQYY | DITLPLLLDS | LQKMQEEMIK | ALKGIFDEYS  | QITSLVTEEI | 240 |
| 241                                     | VNVHKEIQMS | VEQIDPSTEY | NNFIDVHRTT | AAKEQEIEFD | TSLLEENENL  | QANEIMWNNL | 300 |
| 301                                     | TAESLQVMLK | TLAEELMQTQ | QMLLNKEEAV | LELEKRIEES | SETCEKSDI   | VLLLSQKQAL | 360 |
| 361                                     | EELKQSVQQL | RCTEAKFSAQ | KELLEQKVQE | NDGKEPPPVV | NYEEDARSVT  | SMERKERLSK | 420 |
| 421                                     | FESIRHSIAG | IIRSPKSAVG | SSALSDMISI | SEKPLAEQDW | YHGAIPIREA  | QELLKKQGDF | 480 |
| 481                                     | LVRESHGKPG | EYVLSVYSDG | QRRHFIIQYV | DNMYRFEGTG | FSNIPQLIDH  | HYTTKQVITK | 540 |
| 541                                     | KSGVLLNPI  | PKDKKWILSH | EDVILGELLG | KGNFGEVYKG | TLKDKTSVAV  | KTCKEDLPQE | 600 |
| 600                                     | LKIKFLQEA  | ILKQYDHPNI | VKLIGVCTQR | QPVYIIMELV | SGGDFLTLFLR | RKKDELKLNQ | 660 |
| 661                                     | LVKFSLDAAA | GMLYLESKNC | IHRDLAARNC | LVGENNVLKI | SDFGMSRQED  | GGVYSSSGLK | 720 |
| 721                                     | QIPIKWTAPE | ALNYGRYSSE | SDVWSFGILL | WETFSLGVC  | YPGMTNQAR   | EQVERGYRMS | 780 |
| 781                                     | APQHCPEDIS | KIMMKWDYK  | PENRPFSEL  | QKELTIKRK  | LT          |            | 840 |

blue: FER sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](#) accession number NP\_005237.1

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.