

## TIE2 wt

TEK receptor tyrosine kinase

Recombinant Human Active Protein Kinase

HGNC Symbol: TEK

Synonyms: CD202b, TIE-2, VMCM1

Product No.: 0861-0000-1

Lot: 008

**Description:** Human TIE2, C-terminal fragment, amino acids Q<sub>771</sub>-A<sub>1124</sub> (as in [NCBI/Protein](#) entry NP\_000450.1), N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

**Product identity:** TIE2 wt Lot 008, was confirmed as TIE2 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW**<sub>Fusion Protein</sub>: 68,979 Da Da

**Expression host:** Sf9 insect cells

**Purification:** GST-Affinity Chromatography

**Activation:** in vitro auto activation

**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:** 0.143 µg/µl

(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

### Biochemical Parameters:

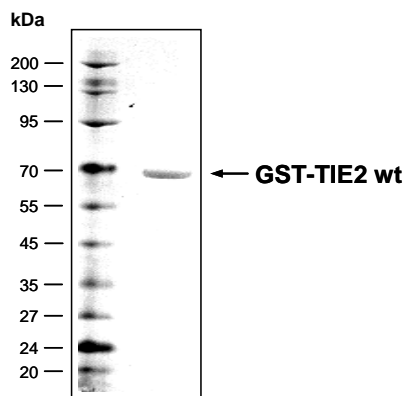
Specific kinase activity (P<sub>i</sub> transfer): 97 pmol/µg × min

ATP-K<sub>M</sub>: 2.2 µM

### Additional assay technology:

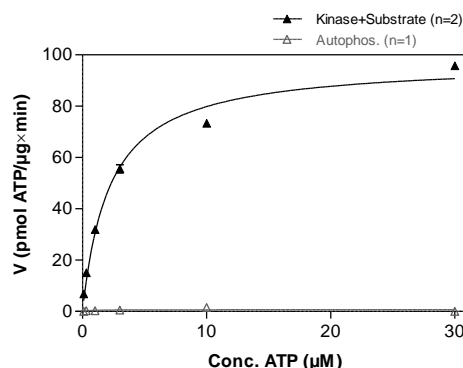
TIE2 wt Lot 008 was also successfully tested by ProQinase for the use with the ADP-Glo™ Kinase assay from Promega  
ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

### TIE2 wt Lot 008: Coomassie stain



2.0 µg GST-TIE2 wt

### TIE2 wt Lot 008: 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: Poly(Ala,Glu,Lys,Tyr)<sub>6:2:5:1</sub> 40 µg/ml
  - Kinase: 1 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

## TIE2

Product No.: 0861-0000-1

GST-TIE2 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDKVLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPOID	KYLKSSKYIA	WPLQGQWQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVLFCG	240
241	PLAMLQLKRA	NVQRMAQAF	QNVREPAVQ	FNSGTLALNR	KVKNNPDPTI	YPVLDWNDIK	300
301	FQDVIGEGNF	GQVLKARIKK	DGLRMDAAIK	RMKEYASKDD	HRDFAGELEV	LCKLGHPNI	360
361	INLLGACEHR	GYLylaIEYA	PHGNLLDFLR	KSRVLETDPA	FAIANSTAST	LSSQQLLHFA	420
421	ADVARGMDYL	SQKQFIHRDL	AARNILVGEN	YVAKIADFGL	SRGQEVYVKK	TMGRLPVRWM	480
481	AIESLNYSVY	TTNSDVWSYG	VLLWEIVSLG	GTPYCGMTCA	ELYEKLPGY	RLEKPLNCDD	540
541	EVYDLMRQCW	REKPYERPSF	AQILVSLNRM	LEERKTYVNT	TLYEKFTYAG	IDCSAEAA	600

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: TIE2 fragment

TIE2 wt <sup>1</sup> Amino Acid Sequence							
1	MDSLASLVLC	GVSLLLSGTV	EGAMDLILIN	SLPLVSDAET	SLTCIASGWR	PHEPITIGRD	60
61	FEALMNQHOD	PLEVTQDVTR	EWAKKVVWKR	EKASKINGAY	FCEGRVRGEA	IRIRTMKMRQ	120
121	QASFLPATLT	MTVDKGDNVN	ISFKKVLIKE	EDAVIYKNGS	FIHSVPRHEV	PDILEVHLP	180
181	AQPQDAGVYS	ARYIGGNLFT	SAFTRLIVRR	CEAQKWGPEC	NHLCTACMNN	GVCHEDTGEC	240
241	ICPPGFMGRT	CEKACELHTF	GRTCKERCSCG	QEGCKSYVFC	LPDPYGCSCA	TGWKGLQCNE	300
301	ACHPGFYGPD	CKLRCSNNG	EMCDRFQGL	CSPGWQGLQC	EREGIPRMTF	KIVDLPDIE	360
361	VNSGKFNPIC	KASGWPLPTN	EEMTLVKPDG	TVLHPKDFNH	TDHFSVAIFT	IHRILPPDSG	420
421	VWVCSVNTVA	GMVEKPFNIS	VKVLPKPLNA	PNVIDTGHN	AVINISSEPY	FGDGPIKSKK	480
481	LLYKPVNHYE	AWQHIQVTNE	IVTLNLYEPR	TEYELCVQLV	RRGEGGEGHP	GPVRRFTTAS	540
541	IGLPPPRGLN	LLPKSQTTLN	LTWQPIFPSS	EDDFYVEVER	RSVQKSDQQN	IKVPGNLTST	600
600	LLNNLHPREQ	YVVRARVNTK	AQGEWSEDLT	AWTLSDILPP	QPENIKISNI	THSSAVISWT	660
661	ILDGYSISSI	TIRYKVQGN	EDQHVDVKIK	NATIIQYQLK	GLEPETAYQV	DIFAENNIGS	720
721	SNPAFSHEL	TLPESQAPAD	LGGGKMLLIA	ILGSAGMTCL	TVLLAFLIIL	QLKRANVQRR	780
781	MAQAFQNVRE	EPAVQFNSGT	LALNRKVKNN	PDPTIYPVLD	WNDIKFQDVI	GEGNFGQVLK	840
841	ARIKDGRLM	DAAIKRMKEY	ASKDDHRDFA	GELEVLCKLG	HHPNIINLLG	ACEHRGYLYL	900
901	AIEYAPHGNL	LDFLRKSRL	ETDPAFAIAN	STASTLSSQQ	LLHFAADVAR	GMDYLSQKQF	960
961	IHRDLAARNI	LVGENYVAKI	ADFGLSRGQE	VYVKKTMGRL	PVRWMAIESL	NYSVYTTNSD	1020
1021	VWSYGVLLWE	IVSLGGTPYC	GMTCAELYEK	LPQGYRLEKP	LNCDDDEVYDL	MRQCWREKPY	1080
1081	ERPSFAQILV	SLNRMLEERK	TYVNTTLYEK	FTYAGIDCSA	EEAA		1140

blue: TIE2 sequence expressed in recombinant protein

<sup>1</sup>NCBI/Protein accession number NP\_000450.1

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.