

## ProQinase™ EGF-R d752-759

epidermal growth factor receptor

Recombinant Human Active Protein Kinase

Synonyms: ERBB, ERBB1

Product No.: 1040-0000-1

Lot: 001

**Description:** Human EGF-R, C-terminal fragment, amino acids H<sub>672</sub>-A<sub>1210</sub> (as in GenBank entry NM\_005228.3), amino acids 752-759 deleted, N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

**Product identity:** EGF-R d752-759 Lot 001, was confirmed as EGF-R by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>Fusion Protein</sub>:** 88,303 Da

**Expression:** Baculovirus infected Sf9 cells

**Purification:** GST-Affinity Chromatography

**Storage buffer:** 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20% glycerol

**Storage temperature:** -80°C  
Avoid repeated freeze-thaw cycles!

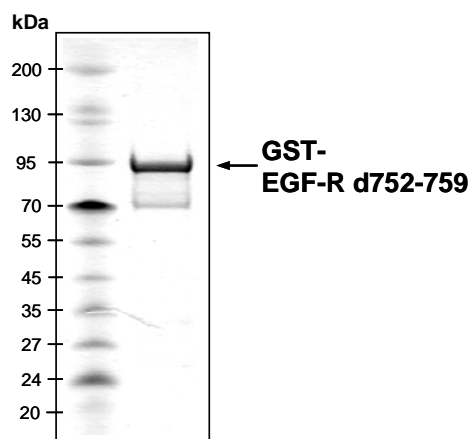
**Protein concentration:** 0.151 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

### 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(Glu:Tyr)<sub>4:1</sub> (Sigma P-0275), 5.0 µg / ml
  - EGF-R d752-759: 1.0 µg / ml
- Filter binding assay
  - MSFC membrane (Millipore)

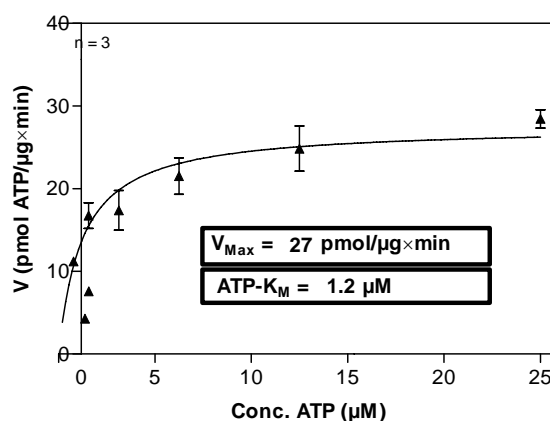
**Specific activity:** 27 pmol/µg×min

### EGF-R d752-759 Lot 001: Coomassie stain



2,0 µg GST-EGF-R d752-759

### EGF-R d752-759 Lot 001: Determination of V<sub>max</sub> and K<sub>m</sub> value for ATP



# ProQinase™ EGF-R d752-759

Product No.: 1040-0000-1

Recombinant Proteins

## EGF-R d752-759 Recombinant Fusion Protein Amino Acid Sequence

1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYE	HLYERDEGDK	WRNKKFELGL	EFNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RDSL <del>EVLFQ</del>	240
241	PLAMGHIVRK	RTLRRLLQER	ELVEPLTPSG	EAPNQALLRI	LKETEFKIK	VLGSGAFGTV	300
301	YKGLWIPEGE	KVKIPVAIKE	LRE <del>ATL</del> DEAY	VMASVDNPHV	CRLLGICLTS	TVQLITQLMP	360
361	FGCLLDYVRE	HKDNIGSQYL	LNWCVQIAKG	MNYLEDRLV	HRDLAARNVL	VKTPQHVKIT	420
421	DFGLAKLLGA	EEKEYHAEGG	KVPIKWMAL	SILHRIYTHQ	SDVWSYGVTV	WELMTFGSKP	480
481	YDGIPASEIS	SILEKGERLP	QPPICTIDVY	MIMVKCWMID	ADSRPKFREL	IIEFSKMARD	540
541	PQRYLVIQGD	ERMHLPSPTD	SNFYRALMDE	EDMDDVVDAD	EYLIPQQGFF	SSPSTSRTP	600
600	LSSLSATSN	STVACIDRNG	LQSCPIKEDS	FLQRYSSDPT	GALTEDSID	TFLVPPEYIN	660
661	QSVPKRPAGS	VQNPVYHNQP	LNPAPSRDPH	YQDPHSTAVG	NPEYLNTPVQ	TCVNFDFSP	720
721	AHWAQKGS HQ	ISLDNPDYQQ	DFFPKAKPN	GIFKGSTAEN	AEYLRVAPQS	SEFIGA	780

1-218: GST    Red: HIS6-tag    Green: 3C    blue: EGF-R fragment    RED: aa flanking the deletion

## EGF-R wt ORF (taken from GenBank entry NM\_005228.3)

1	MRPSGTAGAA	LLALLAALCP	ASRALEEKKV	CQGTSNKL TQ	LGT FEDHFLS	LQRMFN NCEV	60
61	VLGNLEITYV	QRNYDLSFLK	TIQEVAGYVL	IALNTVERIP	LENLQIIRGN	MYYENS YALA	120
121	VLSNYDANKT	GLKELPMRNL	QEILHGAVRF	SNNPALCNVE	SIQWRDIVSS	DFLSNM SMDF	180
181	QNH LGSCQKC	DPSCPNGSCW	GAGEENCQKL	TKIICAQQCS	GRCRGKSPSD	CCHNQCAAGC	240
241	TGPRES DCLV	CRKFRDEATC	KDTCPPMLY	NPTTYQMDVN	PEGKYSFGAT	CVKCKPRNYV	300
301	VTDHGSCVRA	CGADSYEMEE	DGVRKCKKCE	GPCRKVCNGI	GIGEFKDSLS	INATNIKHF	360
361	NCTSISGDLH	ILPVAFRGDS	FTHTPPLDPQ	ELDILKTVKE	ITGFLLIQAW	PENRTDLHAF	420
421	ENLEIIRGRT	KQHGQFSLAV	VSLNITSLGL	RSLKEISDGD	VIISGNKNLC	YANTINWKKL	480
481	FGTSGQKTKI	ISNRGENSCK	ATGQVCHALC	SPEGCWGPEP	RDCVSCRNV	RGRECVDKCN	540
541	LLEGE PREFV	ENSECIQCHP	ECLPQAMNIT	CTGRGPDNCI	QCAHYIDGPH	CVKTC PAVM	600
600	GENNTLVWKY	ADAGHVCHLC	HPNCTYGCTG	PGLEGCP TNG	PKIPSIATGM	VGALLLLLV	660
661	ALGIGLFMR	RHIVRKRTL	RLQERELVE	PLTPSGEAPN	QALLRILKET	EFKKIKVLGS	720
721	GAFGTVYKGL	WIPEGEKVKI	PVAIKELREA	TSPKANKEIL	DEAYVMASVD	NPHVCRLGI	780
781	CLTSTVQLIT	QLMPFGCLLD	YVREHKDNIG	SQYLLNWCVQ	IAGMNYLED	RRLVHRDLAA	840
841	RNVLVKTPQH	VKITDFGLAK	LLGAEKEYEH	AEGGKVPKW	MALESILHRI	YTHQSDVWSY	900
901	GVTWELMTF	GSKPYDGIPA	SEISSILEKG	ERLPQPPICT	IDVYMIMVKC	WMIDADSRPK	960
961	FRELIIIEFSK	MARDPQRYLV	IQGDERMHL	SPTDSNFYRA	LMDEEDMDDV	VDAEYLIPO	1020
1021	QGGFSSPSTS	RTPLLSLSA	TSNNSTVACI	DRNGLQSCPI	KEDSFLQYS	SDPTGALTE	1080
1081	SIDDTFLPVP	EYINQSVPKR	PAGSVQNPVY	HNQPLNPAPS	RDPHYQDPHS	TAVGNPEYLN	1140
1141	TVQPTCVNST	FDSPAHWAQK	GSHQISLDNP	DYQQDFFPKE	AKPNGIFKGS	TAENAEYLRV	1200
1201	APQSSEFIGA						1260

blue: fragment of EGF-R expressed in fusionprotein    RED: 752-759 deletion