

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 H₆₇₂-A₁₂₁₀ (as in GenBank entry NM_005228.3), amino acids 752-759 deleted, N-terminal GST-HIS₆ 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_{Fusion Protein}: 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!

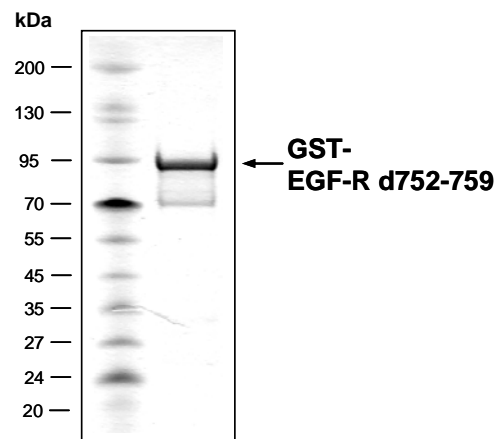
Protein concentration: 0.151 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Determination of K_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: Poly(Glu:Tyr)_{4:1}
(Sigma P-0275), 5.0 µg / ml
 - EGF-R d752-759: 1.0 µg / ml
- Filter binding assay
MSFC membrane (Millipore)

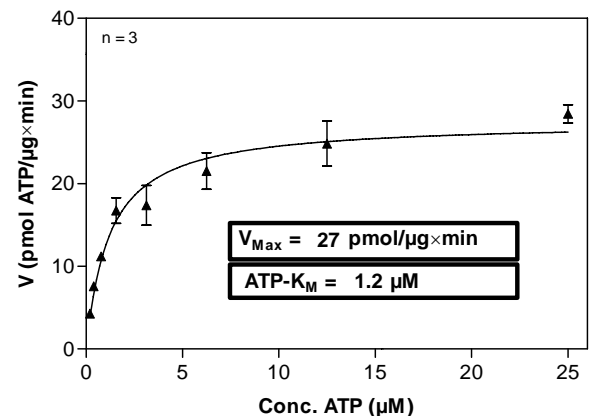
Specific activity: 27 pmol/µg×min

**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_{max} and K_m value for ATP**



This product 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.

EGF-R d752-759

Product No.: 1040-0000-1

EGF-R d752-759 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI	PQID KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHHG	RDSL	240
241	PLAMGHIVRK	RTLRLQLER	ELVEPLTPSG	EAPNQALLRI	LKETEFKKIK	VLGSGAFGTV	300
301	YKGLWIPEGE	KVKIPVAIKE	LREATLDEAY	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	IIEFKSMARD	540
541	PQRYLVIQGD	ERMHLPSPTD	SNFYRALMDE	EDMDDVVDAD	EYLIPQQGFF	SSPSTSRTPL	600
600	LSSLSATSNN	STVACIDRNG	LQSCPIKEDS	FLQRYSSDPT	GALTEDSIDD	TFLPVPEYIN	660
661	QSVPKRPAGS	VQNPVYHNQP	LNPAPSRDPH	YQDPHSTAVG	NPEYLN	TCVNSTFDS	720
721	AHWAQKGS	ISLDNPDYQQ	DDFFPKEAKPN	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	CQGTSNKLTQ	LGTFEDHFLS	LQRMFN	60
61	VLGNLEITYV	QRNYDLSFLK	TIQEVAGYVL	IALNTVERIP	LENLQIIRGN	MYYENS	120
121	VLSNYDANKT	GLKELPMRNL	QEILHGAVRF	SNNPALCNVE	SIQWRDIVSS	DFLSN	180
181	QNHGSCQKC	DPSCPNGSCW	GAGEENCQKL	TKIICAQQCS	GRCRGKSPSD	CCHNQ	240
241	TGPRES	DCLV CRKFRDEATC	KDTCPLMLY	NPTTYQMDVN	PEGKYSFGAT	CVKKC	300
301	VDHGSCVRA	CGADSYEMEE	DGVRKCKKCE	GPCRKVCNGI	GIGEFKDSLS	INATN	360
361	NCTSISGDLH	ILPVAFRGDS	FTHTPPLDPQ	ELDILKTVKE	ITGFLLIQAW	PENRT	420
421	ENLEIIRGRT	KQHGQFSLAV	VSLNITSLGL	RSLKEISDGD	VIISGNKNLC	YANTIN	480
481	FGTSGQKTKI	ISNRGENSCK	ATGQVCHALC	SPEGCWGPEP	RDCVSCRNV	RGREC	540
541	LLEGE	PREFV ENSECIQHP	ECLPQAMNIT	CTGRGPDNCI	QCAHYIDGPH	CVKTC	600
600	GENNTLVWKY	ADAGHVCHLC	HPNCTYGCTG	PGLEGCP	TNG PKIPSIATGM	VGALL	660
661	ALGIGLFMR	RHIVRKRTL	RLLQERELVE	PLTPSGEAPN	QALLRILKET	EFKKIK	720
721	GAFGTVYKGL	WIPEGEKVKI	PVAIKELREA	TSPKANKEIL	DEAYVMASVD	NPHVC	780
781	CLTSTVQLIT	QLMPFGCLLD	YVREHKDNIG	SOYLLNWCVQ	IAKGMNYLED	RRLVHR	840
841	RNVLVKTPQH	VKITDFGLAK	LLGAEKEYH	AEGGKVP	IKW MALESILHRI	YTHQSD	900
901	GVTVWELMTF	GSKPYDGIPA	SEISSILEKG	ERLPQPPICT	IDVYMIMVKC	WMIDADR	960
961	FRELIIEFSK	MARDPQRYLV	IQDERMHL	P SPTDSNFYRA	LMDEEDMDDV	VDADEY	1020
1021	QGFFSSPSTS	RTPLLSSLSA	TSNNSTVACI	DRNGLQSCPI	KEDSFLQRY	SDPTGAL	1080
1081	SIDDTFLPVP	EYINQSVPKR	PAGSVQNPVY	HNQPLNPAPS	RDPHYQDPHS	TAVGNPE	1140
1141	TVQPTCVNST	FDSPA	HWAQK GSHQISLDNP	DYQQDFFPKE	AKPNGIFKGS	TAENAEY	1200
1201	APQSSEFIGA						1260

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

This product 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.