

# EGF-R d747-749/A750P

epidermal growth factor receptor

Recombinant Human Active Protein Kinase

Synonyms: ERBB, ERBB1

Product No.: 1038-0000-1

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 747-749 deleted, A750P mutant, N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

**Product identity:** EGF-R d747-749/A750P Lot 001, was confirmed as EGF-R by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW**<sub>Fusion Protein</sub>: 88,799 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!

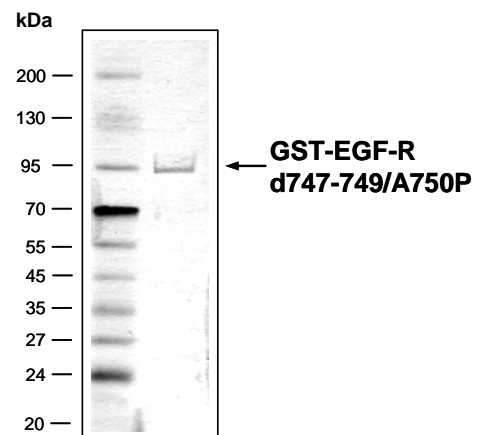
**Protein concentration:** 0.150 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**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 d747-749/A750P: 1.0 µg / ml
- Filter binding assay  
MSFC membrane (Millipore)

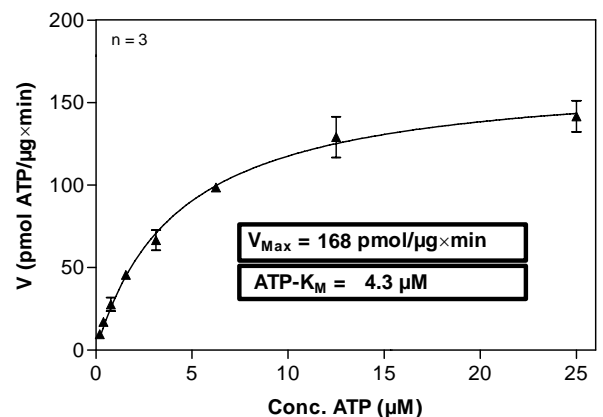
**Specific activity:** 168 pmol/µg×min

**EGF-R d747-749/A750P Lot 001:  
Coomassie stain**



1.0 µg GST-EGF-R d747-749/A750P

**EGF-R d747-749/A750P Lot 001:  
Determination of V<sub>max</sub> and K<sub>m</sub> 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 d747-749/A750P

Product No.: 1038-0000-1

EGF-R d747-749/A750P Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVLFGQ	240
241	PLAMGHIVRK	<u>RTLRLQLQER</u>	<u>ELVEPLTPSG</u>	<u>EAPNQALLRI</u>	<u>LKETEFKKIK</u>	<u>VLGSGAFGTV</u>	300
301	<u>YKGLWIPEGE</u>	<u>KVKIPVAIKE</u>	<u>PTSPKANKEI</u>	<u>LDEAYVMASV</u>	<u>DNPHVCRLLG</u>	<u>ICLTSTVQLI</u>	360
361	<u>TQLMPFGCLL</u>	<u>DYVREHKDNI</u>	<u>GSQYLLNWCV</u>	<u>QIAKGMNYLE</u>	<u>DRRLVHRDLA</u>	<u>ARNVLVKTPQ</u>	420
421	<u>HVKITDFGLA</u>	<u>KLLGAEKEY</u>	<u>HAEGGKVIK</u>	<u>WMALESILHR</u>	<u>IYTHQSDVWS</u>	<u>YGVTVWELMT</u>	480
481	<u>FGSKPYDGIP</u>	<u>ASEISSILEK</u>	<u>GERLPQPPIC</u>	<u>TIDVYMIMVK</u>	<u>CWMIDADSRP</u>	<u>KFRELIIEFS</u>	540
541	<u>KMARDPQRYL</u>	<u>VIQGDERMHL</u>	<u>PSPTDSNFYR</u>	<u>ALMDEEDMDD</u>	<u>VVDADEYLIP</u>	<u>QQGFFSSPST</u>	600
600	<u>SRTPLLSLS</u>	<u>ATSNNSTVAC</u>	<u>IDRNLQSCP</u>	<u>IKEDSFLQRY</u>	<u>SSDPTGALTE</u>	<u>DSIDDTFLPV</u>	660
661	<u>PEYINQSVPK</u>	<u>RPAGSVQNPV</u>	<u>YHNQPLNPAP</u>	<u>SRDPHYQDPH</u>	<u>STAVGNPEYL</u>	<u>NTVQPTCVNS</u>	720
721	<u>TFDSPAHWAQ</u>	<u>KGSHQISLDN</u>	<u>PDYQQDFFPK</u>	<u>EAKPNGIFKG</u>	<u>STAENAEYLR</u>	<u>VAPQSSEFIG</u>	780
781	<u>A</u>						840

1-218: GST **Red:** HIS6-tag **Green:** 3C **blue:** EGF-R fragment **underlined:** aa flanking the deletion **red:** point mutation

EGF-R wt ORF (taken from GenBank entry NM_005228.3)							
1	MRPSGTAGAA	LLALLAALCP	ASRALEEKKV	CQGTSNKLTQ	LGTTFEDHFLS	LQRMFNCEV	60
61	VLGNLEITYV	QRNYDLSFLK	TIQEVAGYVL	IALNTVERIP	LENLQIIRGN	MYYENSYALA	120
121	VLSNYDANKT	GLKELPMRNL	QEILHGAVRF	SNNPALCNVE	SIQWRDIVSS	DFLSNMSMDF	180
181	QNHGSCQKC	DPSPNGSCW	GAGEENCQKL	TKIICAQQCS	GRCRGKSPSD	CCHNQCAAGC	240
241	TGPRESCLV	CRKFRDEATC	KDTCPLMLY	NPTTYQMDVN	PEGKYSFGAT	CVKKCPRNYV	300
301	VDHGSCVRA	CGADSYEMEE	DGVRKCKKCE	GPCRKVCNGI	GIGEFKDSLS	INATNIKHFK	360
361	NCTSISGDLH	ILPVAFRGDS	FTHTPPLDPQ	ELDILKTVKE	ITGFLLIQAW	PENRTDLHAF	420
421	ENLEIIRGRT	KQHGQFSLAV	VSLNITSLGL	RSLKEISDGD	VIISGNKNLC	YANTINWKKL	480
481	FGTSGQTKI	ISNRGENSCK	ATGQVCHALC	SPEGCWPEP	RDCVSCRNVS	RGRECVDKCN	540
541	LLEGEPREFV	ENSECIQCHP	ECLPQAMNIT	CTGRGPDNCI	QCAHYIDGPH	CVKTCPAGVM	600
600	GENNTLVWKY	ADAGHVCHLC	HPNCTYGCTG	PGLEGCPNTG	PKIPSIATGM	VGALLLLLTV	660
661	ALGIGLFMRR	<u>RHIVRKRTL</u>	<u>RLQERELVE</u>	<u>PLTPSGEAPN</u>	<u>QALLRILKET</u>	<u>EFKKIKVLGS</u>	720
721	<u>GAFGTVYKGL</u>	<u>WIPEGEKVKI</u>	<u>PVAIKELREA</u>	<u>TSPKANKEIL</u>	<u>DEAYVMASVD</u>	<u>NPHVCRLGII</u>	780
781	<u>CLTSTVQLIT</u>	<u>QLMPFGCLLD</u>	<u>YVREHKDNIG</u>	<u>SQYLLNWCVQ</u>	<u>IAKGMNYLED</u>	<u>RRLVHRDLAA</u>	840
841	<u>RNVLVKTPQH</u>	<u>VKITDFGLAK</u>	<u>LLGAEKEYH</u>	<u>AEGGKVIKW</u>	<u>MALESILHRI</u>	<u>YTHQSDVWSY</u>	900
901	<u>GVTVWELMTF</u>	<u>GSKPYDGIPA</u>	<u>SEISSILEKG</u>	<u>ERLPQPPICT</u>	<u>IDVYMIMVKC</u>	<u>WMIDADSRPK</u>	960
961	<u>FRELIIEFSK</u>	<u>MARDPQRYLV</u>	<u>IQGDERMHL</u>	<u>SPTDSNFYRA</u>	<u>LMDEEDMDDV</u>	<u>VDADEYLIPQ</u>	1020
1021	<u>QQGFFSSPSTS</u>	<u>RTPLLSLSA</u>	<u>TSNNSTVACI</u>	<u>DRNLQSCPI</u>	<u>KEDSFLQRY</u>	<u>SDPTGALTE</u>	1080
1081	<u>SIDDTFLPV</u>	<u>EYINQSVPKR</u>	<u>PAGSVQNPVY</u>	<u>HNQPLNPAPS</u>	<u>RDPHYQDPHS</u>	<u>TAVGNPEYLN</u>	1140
1141	<u>TVQPTCVNST</u>	<u>FDSPAHWAQK</u>	<u>GSHQISLDNP</u>	<u>DYQQDFFPKE</u>	<u>AKPNGIFKGS</u>	<u>TAENAEYLRV</u>	1200
1201	<u>APQSSEFIGA</u>						1260

**blue:** fragment of EGF-R expressed in fusionprotein **RED:** 747-749 deletion A750 mutation site

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.