

## AXL

AXL receptor tyrosine kinase

Recombinant Human Active Protein Kinase

HGNC Symbol: AXL

Synonyms: EC, UFO

Product No.: 0525-0000-1

Lot: 003

**Description:** Human AXL, C-terminal fragment, amino acids H<sub>464</sub>-A<sub>885</sub> (as in [NCBI/Protein](#) entry NP\_001690.2), N-terminal GST-HIS<sub>6</sub> fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

**Product identity:** AXL, Lot 003, was confirmed as AXL by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW**<sub>Fusion Protein</sub>: 76,549 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:** 0.228 µg/µl

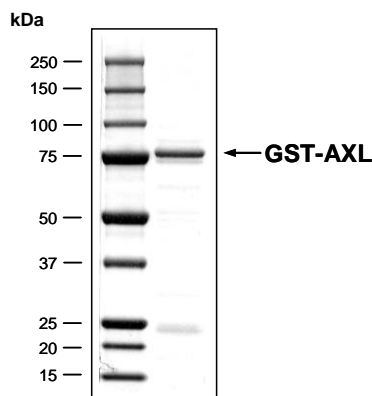
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

### Biochemical Parameters:

Specific kinase activity (P<sub>i</sub> transfer): 26 pmol/µg × min

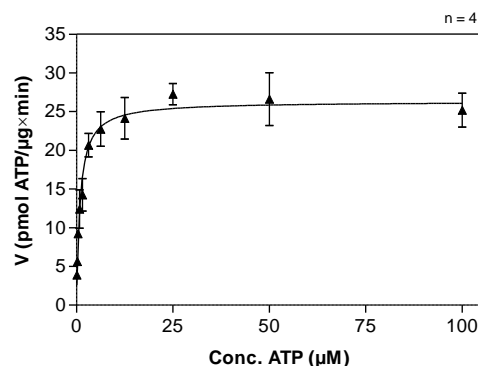
ATP-K<sub>M</sub>: 0.9 µM

### AXL Lot 003: Coomassie stain



2.0 µg GST-AXL

### AXL Lot 003: 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(Glu:Tyr)<sub>4:1</sub>, 2.5 µg/ml
  - Kinase: 1.0 µg/ml
- Filter binding assay
- MSFC membrane (Millipore)

### Additional assay technology:

AXL Lot 003 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.

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.

## AXL

Product No.: 0525-0000-1

GST-AXL Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	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	PMGHHHHHGG	RRRASVAAGI	240
241	LVPRGSPGLD	GICSSHRRRK	ETRYGEVFEP	TVERGELVVR	YRVRKYSRR	TTEATLNSLG	300
301	I SEELKEKLR	DVMVDRHKVA	LGKTLGEGEF	GAVMEGQLNQ	DDSILKVAVK	TMKIAICTRS	360
361	ELEDFLSEAV	CMKEFDHPNV	MRLIGVCFQG	SERESFPAPV	VILPFMKHGD	LHSFLLYSRL	420
421	GDQPVYLP TQ	MLVKFMADIA	SGMEYLSTKR	FIHRDLAARN	CMLNENMSVC	VADFGLSKKI	480
481	YNGDYRQGR	I AKMPVKWIA	IESLADRVYT	SKSDVWSFGV	TMWEIATRQ	TPYPGVENSE	540
541	IYDYLRQGNR	LKQPADCLDG	LYALMSRCWE	LNPQDRPSFT	ELREDLENTL	KALPPAQEPD	600
600	EILYVNMDEG	GGYPEPPGAA	GGADPPTQPD	PKDSCSCLTA	AEVHPAGRYV	LCPSTTPSPA	660
661	QPADRGSPAA	PGQEDGA					720

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

AXL wt <sup>1</sup> Amino Acid Sequence							
1	MAWRCPRMGR	VPLAWCLALC	GWACMAPRGT	QAEESPFVGN	PGNITGARGL	TGTLRCQLQV	60
61	QGEPPEVHWL	RDGQILELAD	STQTQVPLGE	DEQDDWIVVS	QLRITSLQLS	DTGQYQCLVF	120
121	LGHQTFV SQP	GYVGLEGLPY	FLEEPEDRTV	AANTPFNLSC	QAQGPPEPVD	LLWLQDAVPL	180
181	ATAPGHG PQR	SLHVPGLNKT	SSFSCAEHNA	KGVTTSR TAT	ITVLPQQPRN	LHLVSRQPT E	240
241	LEVAWTPGLS	GIYPLTHCTL	QAVLSDDGMG	IQAGEPDPPE	EPLTSQASVP	PHQLRLGSLH	300
301	PHTPYHIRVA	CTSSQG PSSW	THWLPVETPE	GVPLGPPENI	SATRNGSQAF	VHWQEPRAPL	360
361	QGTLLGYRLA	YQGQDTPEVL	MDIGLRQEV T	LELQGDG SVS	NLTVCVAA YT	AAGDGPWSLP	420
421	VPLEAWRPVK	EPSTPAFSWP	WWYVLLGAVV	AAACVLILAL	FLVHRRK KET	RYGEVF EPTV	480
481	ERGELVVRYR	VRKYSRRRT	EATLNSLGIS	EELKEKLRDV	MVDRHKVALG	KTLGEGEF GA	540
541	VMEGQLNQDD	SILKVAVKTM	KIAICTRSEL	EDFLSEAVCM	KEFDHPNVMR	LIGVCFQ GSE	600
600	RESFPAPVVI	LPFMKHGD LH	SFLLYSRLGD	QPVYLP TQML	VKFMADIASG	MEYLSTKR FI	660
661	HRDLAARNCM	LNENMSVCVA	DFGLSKKIYN	GDYRQGR IA	KMPVKWIAIE	SLADRVYTSK	720
721	SDVWSFGVTM	WEIATRQ T P	YPGVENSEIY	DYLRQGNRLK	QPADCLDGLY	ALMSRCWELN	780
781	PQDRPSFTEL	REDLENTLKA	LPPAQEPDEI	LYVNMDEGGG	YPEPPGAAGG	ADPPTQDPK	840
841	DSCSCLTAAE	VHPAGRYVLC	PSTTPSPAQP	ADRGSPAAPG	QEDGA		900

**blue:** AXL sequence expressed in recombinant protein

<sup>1</sup>[NCBI/Protein](#) accession number NP\_001690.2

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.