

## ProQinase™ 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

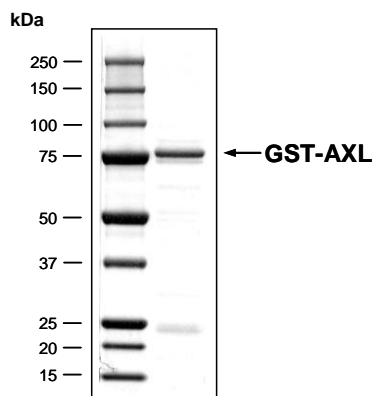
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

### Biochemical Parameters:

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

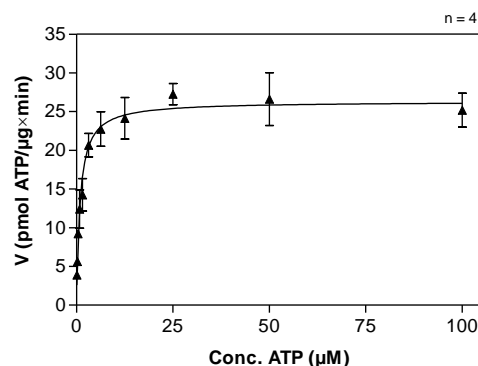
ATP-K<sub>M</sub>: 0.9 µM

### 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 Reaction Biology 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

## ProQinase™ AXL

Product No.: 0525-0000-1

GST-AXL Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSM	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQG WQATF	GGGDHPPKSD	PMG <b>HHHHHG</b>	RRRASVAAGI	240
241	<b>LVPRGS</b> PGLD	GICSSHRRRK	<b>ETRYGEVFEP</b>	<b>TVERGELVVR</b>	<b>YRVRKYSRR</b>	<b>TTEATLNSLG</b>	300
301	<b>I SEELKEKLR</b>	<b>DVMVDRHKVA</b>	<b>LGKTLGEGEF</b>	<b>GAVMEGQLNQ</b>	<b>DDSILKVAVK</b>	<b>TMKIAICTRS</b>	360
361	<b>ELEDFLSEAV</b>	<b>CMKEFDHPNV</b>	<b>MRLIGVCFQG</b>	<b>SERESFPAPV</b>	<b>VILPFMKHGD</b>	<b>LHSFLLYSRL</b>	420
421	<b>GDQP VYLPTQ</b>	<b>MLVKFMADIA</b>	<b>SGMEYLSTKR</b>	<b>FIHRDLAARN</b>	<b>CMLNENMSVC</b>	<b>VADFGLSKKI</b>	480
481	<b>YNGDYRQGR</b>	<b>I AKMPVKWIA</b>	<b>IESLADRVYT</b>	<b>SKSDVWSFGV</b>	<b>TMWEIATRQ</b>	<b>TPYPGVENSE</b>	540
541	<b>IYDYLRQGNR</b>	<b>LKQPADCLDG</b>	<b>LYALMSRCWE</b>	<b>LNPQDRPSFT</b>	<b>ELREDLENTL</b>	<b>KALPPAQEPD</b>	600
600	<b>EILYVNMDEG</b>	<b>GGYPEPPGAA</b>	<b>GGADPPTQPD</b>	<b>PKDSCSCLTA</b>	<b>AEVHPAGRYV</b>	<b>LCPSTTPSPA</b>	660
661	<b>QPADRGSPA</b>	<b>PGQEDGA</b>					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	KGVTTSRTAT	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 Y T	AAGDGPWSLP	420
421	VPLEAWRPVK	EPSTPAFSWP	WWYVLLGAVV	AAACVLILAL	FLV <b>HRRKKET</b>	<b>RYGEVF EPTV</b>	480
481	<b>ERGELVVRYR</b>	<b>VRKYSRRRT</b>	<b>EATLNSLGIS</b>	<b>EELKEKLRDV</b>	<b>MVDRHKVALG</b>	<b>KTLGEGEFGA</b>	540
541	<b>VMEGQLNQDD</b>	<b>SILKVAVKTM</b>	<b>KIAICTRSEL</b>	<b>EDFLSEAVCM</b>	<b>KEFDHPNVMR</b>	<b>LIGVCFQ GSE</b>	600
600	<b>RESFPAPVVI</b>	<b>LPFMKHGD LH</b>	<b>SFLLYSRLGD</b>	<b>QPVYLPTQML</b>	<b>VKFMADIASG</b>	<b>MEYLSTKRFI</b>	660
661	<b>HRDLAARNCM</b>	<b>LNENMSVCVA</b>	<b>DFGLSKKIYN</b>	<b>GDYRQGR IA</b>	<b>KMPVKWIAIE</b>	<b>SLADRVYTSK</b>	720
721	<b>SDVWSFGVTM</b>	<b>WEIATRQTP</b>	<b>YPGVENSEIY</b>	<b>DYLRQGNRLK</b>	<b>QPADCLDGLY</b>	<b>ALMSRCWELN</b>	780
781	<b>PQDRPSFTEL</b>	<b>REDLENTLKA</b>	<b>LPPAQEPDEI</b>	<b>LYVNMDEGGG</b>	<b>YPEPPGAAGG</b>	<b>ADPPTQDPK</b>	840
841	<b>DSCSCLTAAE</b>	<b>VHPAGRYVLC</b>	<b>PSTTPSPAQP</b>	<b>ADRGSPAAPG</b>	<b>QEDGA</b>		900

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

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