

AXL (AXL receptor tyrosine kinase)

Synonyms: EC, UFO

Product No.: 0525-0000-1

Lot: 003

Description: Human AXL, C-terminal fragment, amino acids H₄₆₄-A₈₈₅ (as in GenBank entry NM_001699.3)*, N-terminally fused to GST-HIS₆-Thrombin cleavage site, expressed in Sf9 insect cells

*Sequence may contain documented polymorphisms
Detailed aa-sequence on request

Product identity: AXL, Lot 003, was confirmed as AXL by mass spectroscopy LC-ESI-MS/MS (Protagen AG, Germany)

Theoretical MW_{Fusion Protein}: 76,549 Da

Expression: Baculovirus infected Sf9 cells

Purification: GST-Affinity Chromatography

Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 4 mM reduced glutathione, 20% glycerol

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

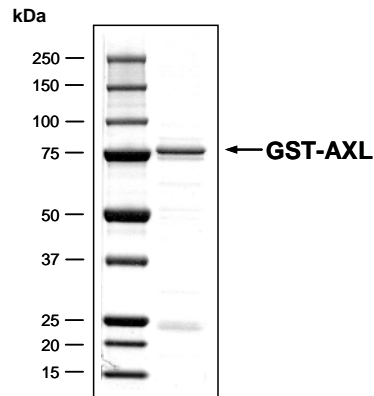
Protein concentration: 0.228 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Method for 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
 - 2.5 µg / 50 µl PEG_{20,000}
 - ATP (variable)
 - Substrate: Poly(Glu:Tyr)_{4:1} (Sigma P-0275), 0.25 µg / 50 µl
 - AXL: 200 ng / 50 µl
- Filter binding assay
MSFC membrane (Millipore)

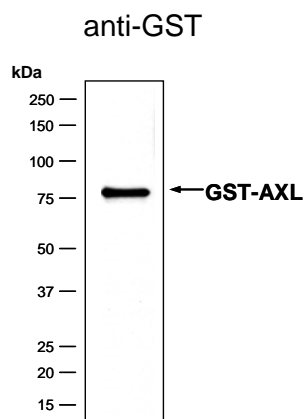
Specific activity: 26 pmol/µg×min

Coomassie stain:



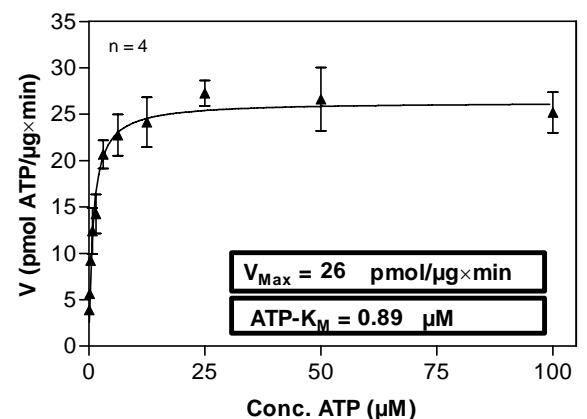
2.0 µg GST-AXL

Western blot analysis:



500 ng GST-AXL

Determination of 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.