

Product datasheet

Human Axl ELISA Kit ab208987

Recombinant SimpleStep ELISA

[6 Images](#)

Overview

Product name Human Axl ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Serum	5			2.3%

Inter-assay

Sample	n	Mean	SD	CV%
Serum	3			2.9%

Sample type Cell culture supernatant, Saliva, Milk, Urine, Serum, Hep Plasma, EDTA Plasma, Cit plasma

Assay type Sandwich (quantitative)

Sensitivity 11 pg/ml

Range 15.6 pg/ml - 1000 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Milk	90	87% - 95%
Urine	108	102% - 115%
Serum	110	109% - 111%
Cell culture media	99	97% - 100%
Hep Plasma	105	105% - 106%
EDTA Plasma	108	107% - 109%

Sample type	Average %	Range
Cit plasma	107	105% - 110%

Assay time

1h 30m

Assay duration

One step assay

Species reactivity

Reacts with: Human

Does not react with: Goat, Cow, Pig

Product overview

Human Axl ELISA Kit (ab208987) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Axl protein in cell culture supernatant, cit plasma, edta plasma, hep plasma, milk, saliva, serum, and urine. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human Axl with 11 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate ([ab203359](#)) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

Notes

Axl is a transmembrane receptor tyrosine kinase on the plasma membrane. The growth factor GAS6 is a ligand for Axl and regulates many physiological processes including cell survival, cell proliferation, migration and differentiation. The antibodies used in this kit were raised against a portion of the extracellular domain of Axl.

Platform

Microplate (12 x 8 well strips)

Properties

Storage instructions

Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Human Axl Capture Antibody	1 x 600µl
10X Human Axl Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent CPI - HAMA Blocker (ab193969)	1 x 6ml

Components	1 x 96 tests
Human Axl Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

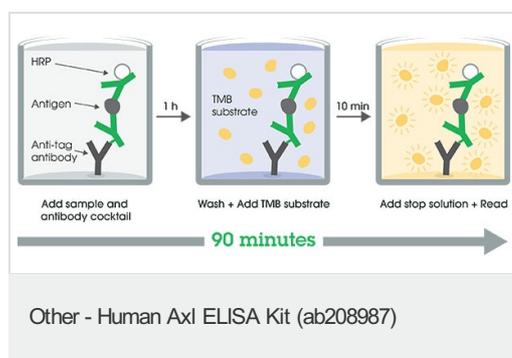
Function May function as a signal transducer between specific cell types of mesodermal origin. In case of filovirus infection, seems to function as a cell entry factor.

Tissue specificity Highly expressed in metastatic colon tumors. Expressed in primary colon tumors. Weakly expressed in normal colon tissue.

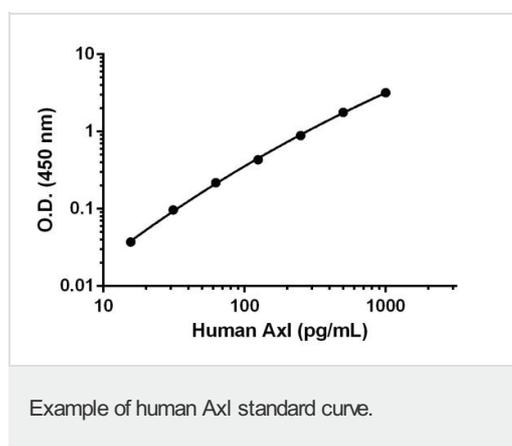
Sequence similarities Belongs to the protein kinase superfamily. Tyr protein kinase family. AXL/UFO subfamily. Contains 2 fibronectin type-III domains. Contains 2 Ig-like C2-type (immunoglobulin-like) domains. Contains 1 protein kinase domain.

Cellular localization Membrane.

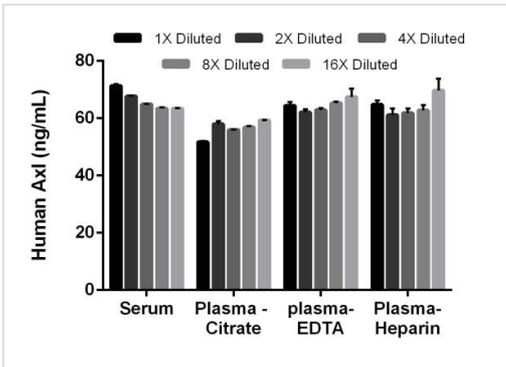
Images



SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.

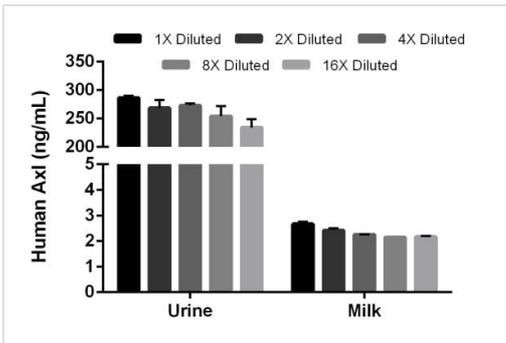


Background-subtracted data values (mean +/- SD) are graphed.



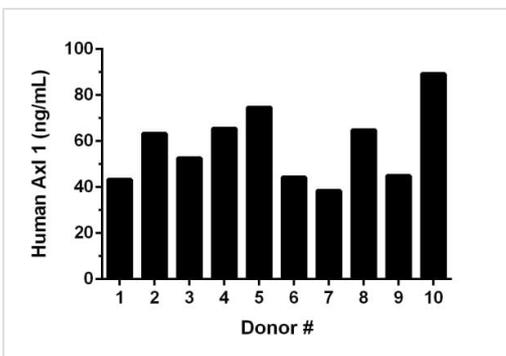
Interpolated concentrations of native Axl in human serum, and plasma samples.

The concentrations of Axl were measured in duplicates, interpolated from the Axl standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 1%, plasma (citrate) 2%, plasma (EDTA) 0.5% and plasma (heparin) 1%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Axl concentration was determined to be 66 ng/mL in serum, 56 ng/mL in plasma (citrate) and 66 ng/mL in plasma (EDTA) and 64 ng/mL in plasma (heparin).



Interpolated concentrations of native Axl in human urine and milk samples.

The concentrations of Axl were measured in duplicates, interpolated from the Axl standard curves and corrected for sample dilution. Undiluted samples are as follows: urine 0.25% and milk 10%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Axl concentration was determined to be 271 ng/mL in urine and 2.38 ng/mL in milk.



Serum from ten individual healthy human female donors was measured in duplicate.

Interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Axl concentration was determined to be 58.1 ng/mL with a range of 37.8 – 89.9 ng/mL.

Powered by
recombinant antibodies



Research with confidence
Consistent and reproducible results



Long-term and scalable supply
Recombinant technology



Success from the first experiment
Confirmed specificity



Ethical standards compliant
Animal-free production

Sandwich ELISA - Human Axl ELISA Kit (ab208987)

To learn more about the advantages of recombinant antibodies see [here](#).

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors