

## Product datasheet

# Human Lipoprotein A ELISA Kit ab108878

[1 References](#) [2 Images](#)

### Overview

**Product name** Human Lipoprotein A ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
--------	---	------	----	-----

Overall 4.6%

Inter-assay

Sample	n	Mean	SD	CV%
--------	---	------	----	-----

Overall 9.5%

**Sample type** Cell culture supernatant, Milk, Urine, Serum, Plasma

**Assay type** Sandwich (quantitative)

**Sensitivity** = 0.7 ng/ml

**Range** 0.78 ng/ml - 50 ng/ml

**Recovery** = 98 %

**Assay time** 4h 00m

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Product overview**

Lipoprotein A Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Lipoprotein A concentrations in plasma, serum, urine, milk, cerebrospinal fluid (CSF) and cell culture supernatants.

A Lipoprotein A specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Lipoprotein A specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the

amount of Lipoprotein A captured in plate.

**The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.**

**Platform** Microplate

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

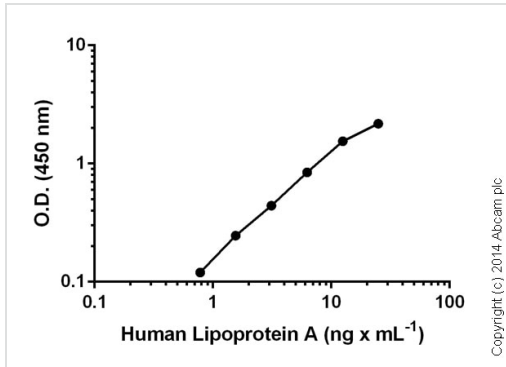
Components	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent M Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human Lipoprotein A Antibody	1 x 120µl
Chromogen Substrate	1 x 8ml
Lipoprotein A Microplate (12 x 8 well strips)	1 unit
Lipoprotein A Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 12ml

**Function** Apo(a) is the main constituent of lipoprotein(a) (Lp(a)). It has serine proteinase activity and is able of autoprolysis. Inhibits tissue-type plasminogen activator 1. Lp(a) may be a ligand for megalin/Gp 330.

**Sequence similarities** Belongs to the peptidase S1 family. Plasminogen subfamily.  
Contains 38 kringle domains.  
Contains 1 peptidase S1 domain.

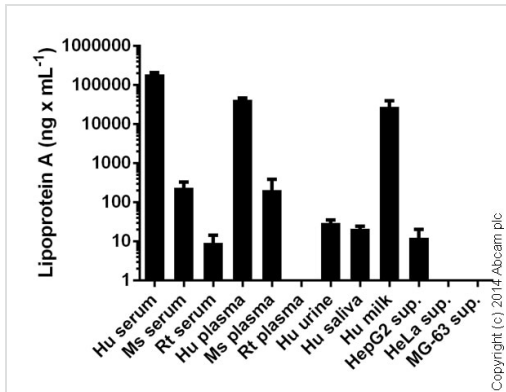
**Post-translational modifications** N- and O-glycosylated. The N-glycans are complex biantennary structures present in either a mono- or disialylated state. The O-glycans are mostly (80%) represented by the monosialylated core type I structure, NeuNAc $\alpha$ 2-3Gal $\beta$ 1-3GalNAc, with smaller amounts of disialylated and non-sialylated O-glycans also detected.

## Images



Standard curve: mean of duplicates (+/- SD) with background reads subtracted

Sandwich ELISA - Lipoprotein A Human ELISA Kit (ab108878)



Lipoprotein A measured in biological fluids and cell culture supernatants showing quantity (ng) per mL of tested sample. Human serum, plasma and milk samples diluted 4000-32000 fold. Human urine and saliva diluted 2-8 fold. Mouse and rat samples diluted 10-100 fold. Cell supernatants diluted 1-10 fold.

Sandwich ELISA - Lipoprotein A Human ELISA Kit (ab108878)

**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