



Product datasheet

Human Lipoprotein A ELISA Kit ab108878

★★★★★ 1 Abreviews 2 References 画像数 2

製品の概要

製品名	Human Lipoprotein A ELISA Kit			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				4%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				9.9%
サンプルの種類	Milk, Urine, Serum, Plasma, Cerebral Spinal Fluid			
アッセイタイプ	Sandwich (quantitative)			
検出感度	= 0.31 ng/ml			
検出範囲	1.563 ng/ml - 50 ng/ml			
添加回収試験	= 102 %			
全工程の試験時間	4h 00m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	Lipoprotein A Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Lipoprotein A concentrations in plasma, serum, urine, milk, CSF, cell culture supernatants, cell lysate, and tissue samples.			

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.

試験プラットフォーム

Microplate

製品の特性

保存方法

Store at -20°C. Please refer to protocols.

内容	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 100µl
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

機能

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

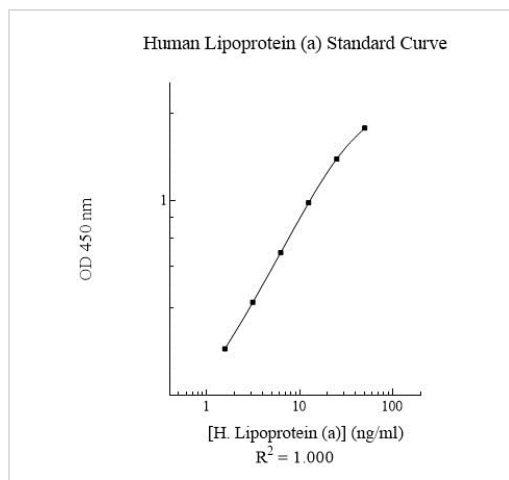
配列類似性

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

翻訳後修飾

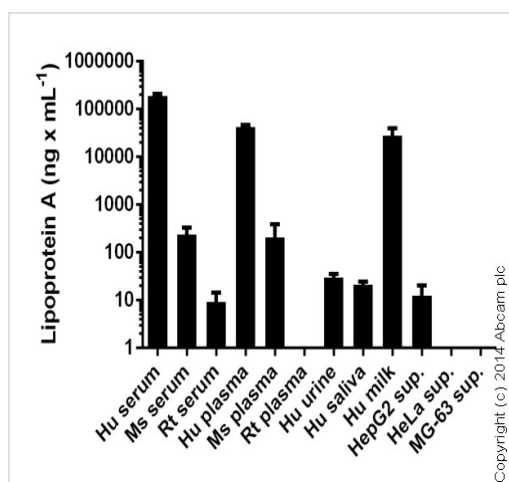
N- and O-glycosylated. The N-glycans are complex biantennary structures present in either a mono- or disialylated state. The O-glycans are mostly (80%) represented by the monosialylated core type I structure, NeuNAcα2-3Galβ1-3GalNAc, with smaller amounts of disialylated and non-sialylated O-glycans also detected.

画像



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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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