

Human Apolipoprotein A-IV ELISA Kit ab214567

リコンビナント SimpleStep ELISA

★★★★★ 1 Abreviews 1 References 画像数 9

製品の概要

製品名	Human Apolipoprotein A-IV 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
	Serum	8		CV%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Serum	3		CV%
サンプルの種類	Cell culture supernatant, Saliva, Urine, Serum, Tissue Extracts, Hep Plasma, EDTA Plasma, Cit plasma			
アッセイタイプ	Sandwich (quantitative)			
検出感度	5.4 pg/ml			
検出範囲	250 pg/ml - 16000 pg/ml			
添加回収試験	特定サンプルでの回収試験			
	サンプルの種類	平均 %	測定範囲	
	Saliva	78	76% - 79%	
	Urine	97	95% - 98%	
	Serum	98	90% - 101%	
	Tissue Extracts	103	102% - 105%	
	Cell culture media	103	102% - 103%	
	Hep Plasma	98	95% - 101%	

サンプルの種類	平均 %	測定範囲
EDTA Plasma	104	102% - 107%
Cit plasma	91	90% - 94%

全工程の試験時間

1h 30m

ステップ

One step assay

種交差性

交差種: Human

非交差種: Cow

製品の概要

Human Apolipoprotein A-IV ELISA Kit (ab214567) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Apolipoprotein A-IV protein in cit plasma, edta plasma, hep plasma, saliva, serum, tissue extracts, urine, and cell culture supernatant. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human Apolipoprotein A-IV with 5.4 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.

特記事項

Human Apolipoprotein A-IV is a 45kDa glycoprotein with activities involving lipid transport. Specifically, Apolipoprotein A-IV is a component of high density lipoprotein (HDL) and chylomicrons. Its primarily synthesized is the intestine in association with lymph chylomicron particles. Although its precise function is not known, Apolipoprotein A-IV is a potent activator of lecithin-cholesterol acyltransferase (LCAT) *in vitro*. Human Apolipoprotein A-IV consists of a 20 amino acid (aa) signal peptide and 376 aa mature chain. Human Apolipoprotein A-IV is 62% identical to both mouse and rat Apolipoprotein A-IV.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

試験プラットフォーム

Pre-coated microplate (12 x 8 well strips)

製品の特性

保存方法

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

内容	1 x 96 tests
10X Wash Buffer PT (ab206977)	1 x 20ml
50X Cell Extraction Enhancer Solution (ab193971)	1 x 1ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml
Antibody Diluent 5BI	1 x 6ml
10X Human Apolipoprotein A-IV Capture Antibody	1 x 600µl
10X Human Apolipoprotein A-IV Detector Antibody	1 x 600µl
Human Apolipoprotein A-IV 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

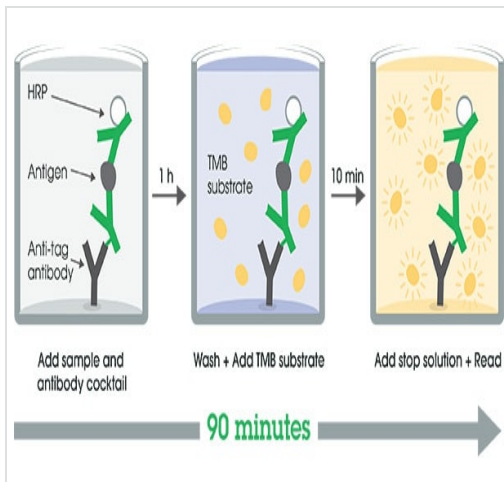
関連性

APOA4 (apolipoprotein A-IV) is a component of HDL and chylomicrons. Its primary site of synthesis is the intestine, in association with lymph chylomicron particles. Although its precise function is not known, APOA4 is a potent activator of lecithin-cholesterol acyltransferase (LCAT) in vitro. In rodents, Apo A-IV inhibits gastric emptying and serves as a satiety factor whose synthesis and secretion are increased by the ingestion of dietary fat. It also possesses anti-inflammatory and antiatherogenic properties

細胞内局在

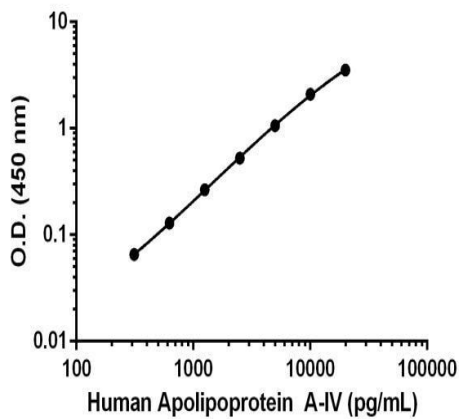
Secreted

画像



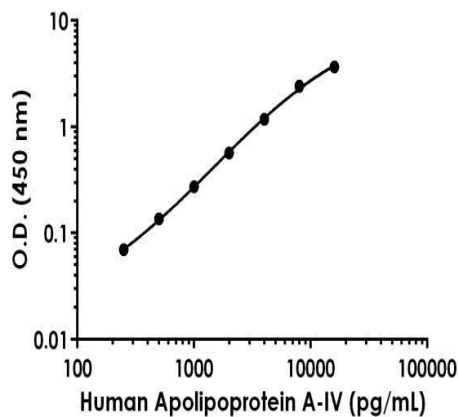
SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.

Other - Human Apolipoprotein A-IV ELISA Kit
(ab214567)



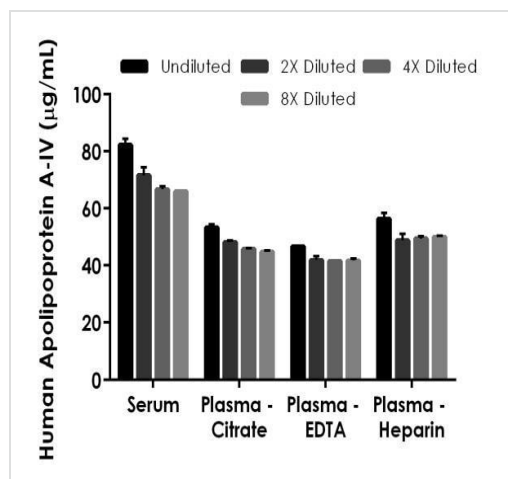
Example of human Apolipoprotein A-IV standard curve in Sample Diluent NS.

Background-subtracted data values (mean \pm SD) are graphed.



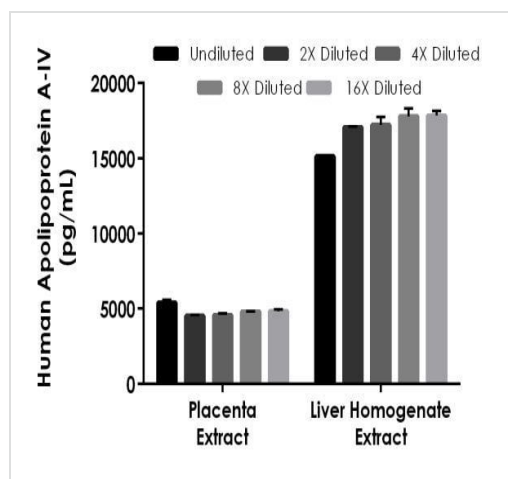
Example of human Apolipoprotein A-IV standard curve in Sample Diluent 1X Cell Extraction Buffer PTR.

Background-subtracted data values (mean \pm SD) are graphed.



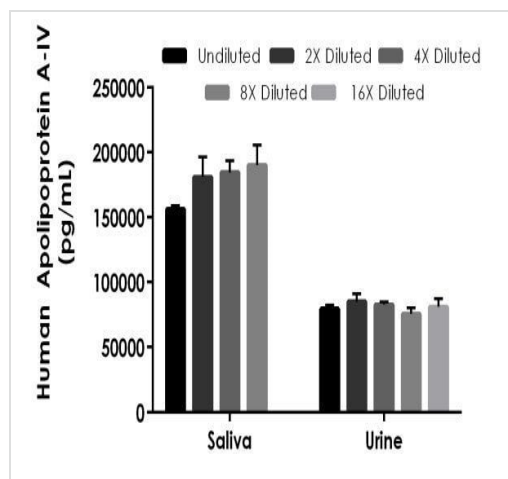
Interpolated concentrations of native Apolipoprotein A-IV in human serum and plasma samples.

The concentrations of Apolipoprotein A-IV were measured in duplicates, interpolated from the Apolipoprotein A-IV standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 1:32,000, plasma (citrate) 1:16,000, plasma (EDTA) 1:16,000, and plasma (heparin) 1:16000. The interpolated dilution factor corrected values are plotted (mean \pm SD, n=2). The mean Apolipoprotein A-IV concentration was determined to be 71.65 μ g/mL in neat serum, 48.04 μ g/mL in neat plasma (citrate), 42.94 μ g/mL in neat plasma (EDTA) and 51.18 μ g/mL in neat plasma (heparin).



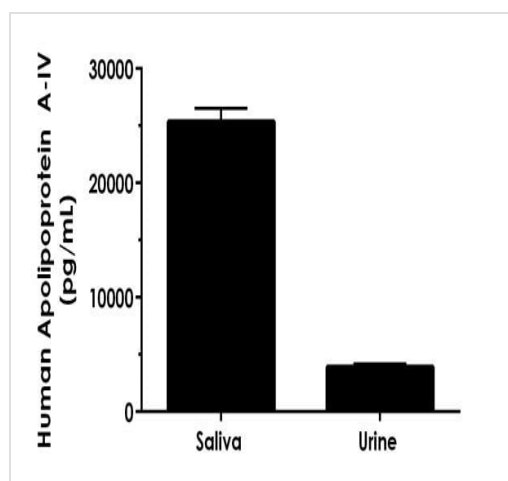
Interpolated concentrations of native Apolipoprotein A-IV in human placenta and human liver homogenate tissue extracts based on 500 μ g/mL and 400 μ g/mL extract loads, respectively.

The concentrations of Apolipoprotein A-IV were measured in duplicate and interpolated from the Apolipoprotein A-IV standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean \pm SD, n=2). The mean Apolipoprotein A-IV concentration was determined to be 4,829 pg/mL in placenta extract and 17,016 pg/mL in liver homogenate extract.



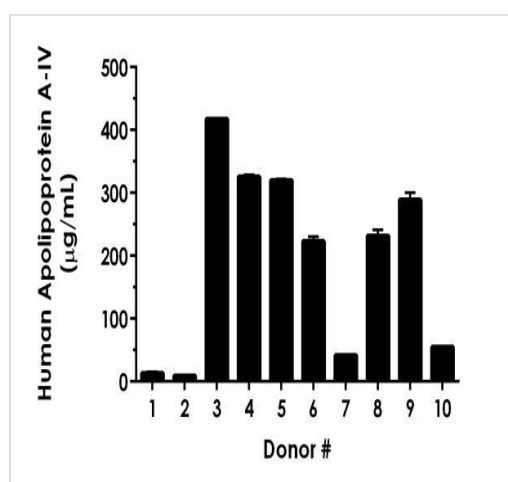
Interpolated concentrations of spiked Apolipoprotein A-IV in human saliva and urine samples.

The concentrations of Apolipoprotein A-IV were measured in duplicates, interpolated from the Apolipoprotein A-IV standard curves and corrected for sample dilution. Undiluted samples are as follows: saliva 6.25% and urine 12.5%. The interpolated dilution factor corrected values are plotted (mean \pm SD, $n=2$). The mean Apolipoprotein A-IV concentration was determined to be 177,941 pg/mL in neat saliva (spiked) and 80,680 pg/mL in neat urine (spiked).



Interpolated concentrations of native Apolipoprotein A-IV in human saliva and urine samples.

Saliva was analyzed at 6.25% and urine at 12.5%. The concentrations of Apolipoprotein A-IV were measured in duplicates, interpolated from the Apolipoprotein A-IV standard curves and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean \pm SD, $n=2$). The mean Apolipoprotein A-IV concentration was determined to be 25,370 pg/mL in neat saliva (native) and 3903 pg/mL in neat urine (native).



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

Interpolated dilution factor corrected values are plotted (mean \pm SD, $n=2$). The mean human Apolipoprotein A-IV concentration was determined to be 192.8 μ g/mL with a range of 9.2 – 418.0 μ g/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 Apolipoprotein A-IV
ELISA Kit (ab214567)

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.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