

Human Apolipoprotein H ELISA Kit (APOH) αb108814

2 Images

Overview

Product name	Human Apolipoprotein H ELISA Kit (APOH)			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			CV%
				3.5%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			CV%
				9.7%
Sample type	Saliva, Milk, Urine, Serum, Plasma			
Assay type	Sandwich (quantitative)			
Sensitivity	= 0.13 ng/ml			
Range	0.313 ng/ml - 20 ng/ml			
Recovery	98 %			
Assay time	4h 00m			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	Abcam's Apolipoprotein H (APOH) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Apolipoprotein H concentrations in Human plasma, serum, saliva, urine, milk and cell culture samples.			

An Apolipoprotein H specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an Apolipoprotein H specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Complex 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 Apolipoprotein H 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

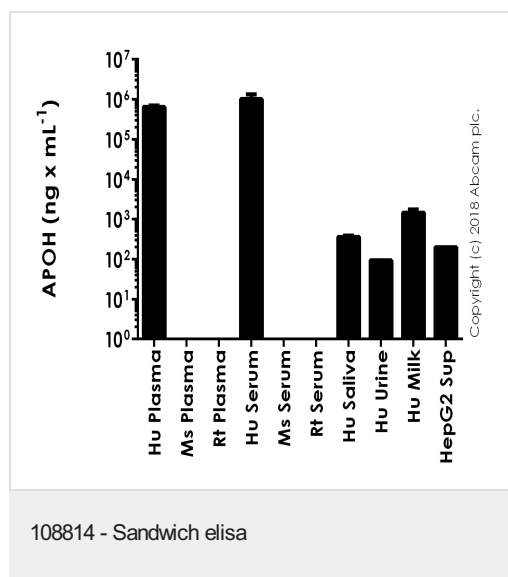
Properties

Storage instructions

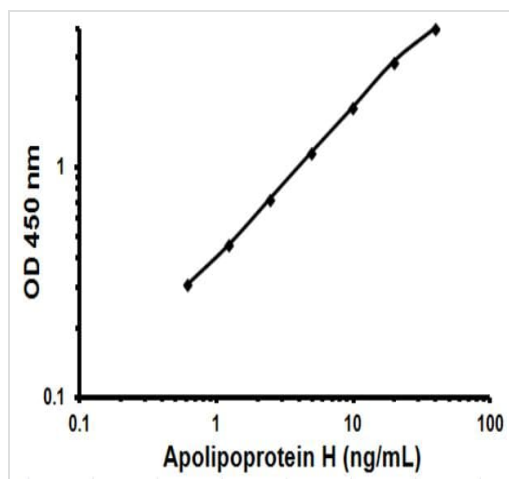
Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human Apolipoprotein H Antibody	1 x 120µl
Apolipoprotein H Microplate (12 x 8 well strips)	1 unit
Apolipoprotein H Standard (Lyophilized)	1 x 24ng
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml

Images



Apolipoprotein H measured in biological fluids with background signal subtracted (duplicates +/- SD). The samples were tested in a range of dilutions (human serum and plasma at 1:10000-1:200000; milk at 1:100-1:1000 and the rest 1:1-1:50).



Representative Standard Curve using ab108814

Typical Standard Curve

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