

Human Apolipoprotein H ELISA Kit (APOH) ab108814

2 图像

概述

产品名称	人Apolipoprotein H ELISA试剂盒(APOH)																				
检测方法	Colorimetric																				
精确度	<div>批次内</div> <table><tr><th>样品</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>Overall</td><td></td><td></td><td></td><td>3.5%</td></tr></table> <div>批次间</div> <table><tr><th>样品</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>Overall</td><td></td><td></td><td></td><td>9.7%</td></tr></table>	样品	n	Mean	SD	CV%	Overall				3.5%	样品	n	Mean	SD	CV%	Overall				9.7%
样品	n	Mean	SD	CV%																	
Overall				3.5%																	
样品	n	Mean	SD	CV%																	
Overall				9.7%																	
样品类型	Saliva, Milk, Urine, Serum, Plasma																				
检测类型	Sandwich (quantitative)																				
灵敏度	= 0.13 ng/ml																				
范围	0.313 ng/ml - 20 ng/ml																				
回收率	98 %																				
检测时间	4h 00m																				
实验步骤	Multiple steps standard assay																				
种属反应性	与反应: Human																				
产品概述	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.

平台

Microplate

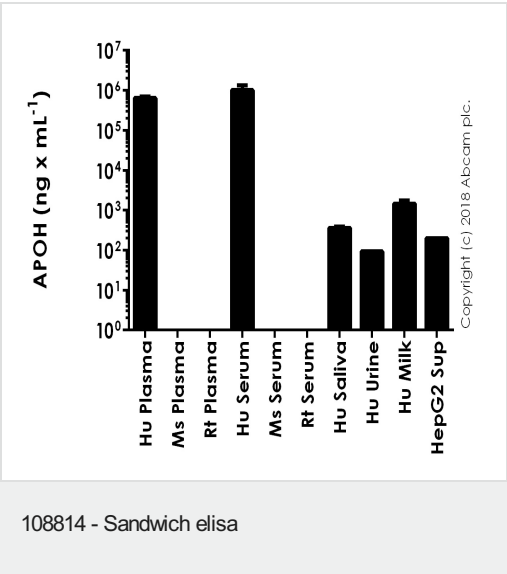
性能

存放说明

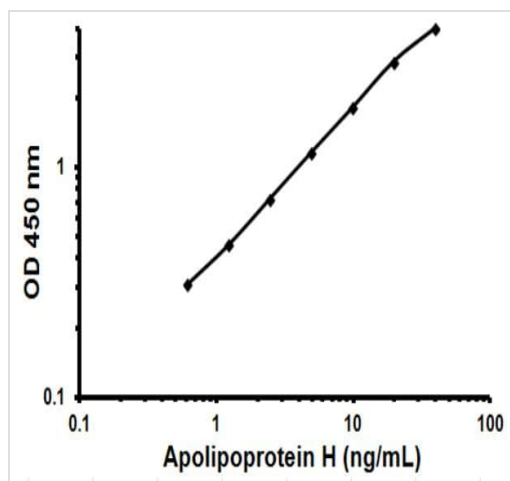
Store at -20°C. Please refer to protocols.

组件	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

图片



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.cn/abpromise> or contact our technical team.

### Terms and conditions

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