

Human Apolipoprotein B ELISA Kit (APOB) ab108807

★★★★★ 1 Abreviews 11 References 2 图像

概述

产品名称 人Apolipoprotein B ELISA试剂盒(APOB)

检测方法 Colorimetric

批次内				
样品	n	Mean	SD	CV%
Overall				6.4%

批次间				
样品	n	Mean	SD	CV%
Overall				10.4%

样品类型 Cell culture supernatant, Serum, Plasma, Cerebral Spinal Fluid

检测类型 Sandwich (quantitative)

灵敏度 = 1.7 ng/ml

范围 3.906 ng/ml - 250 ng/ml

回收率 97 %

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 B (APOB) Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Apolipoprotein B concentrations in Human plasma, serum, CSF and cell culture samples.

An Apolipoprotein B specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an Apolipoprotein B 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 B captured in plate.

Get results in 90 minutes with Human Apolipoprotein B ELISA Kit ([ab190806](#)) from our SimpleStep ELISA® range.

**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 M Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human Apolipoprotein B Antibody	1 x 120µl
Apolipoprotein B Microplate (12 x 8 well strips)	1 unit
Apolipoprotein B Standard	2 vials
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml

功能 Apolipoprotein B is a major protein constituent of chylomicrons (apo B-48), LDL (apo B-100) and VLDL (apo B-100). Apo B-100 functions as a recognition signal for the cellular binding and internalization of LDL particles by the apoB/E receptor.

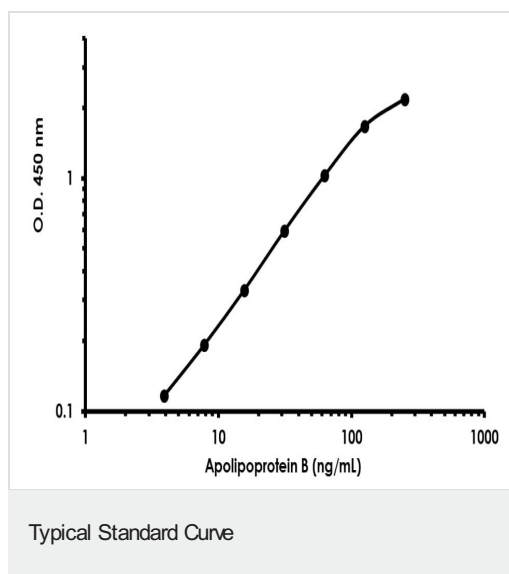
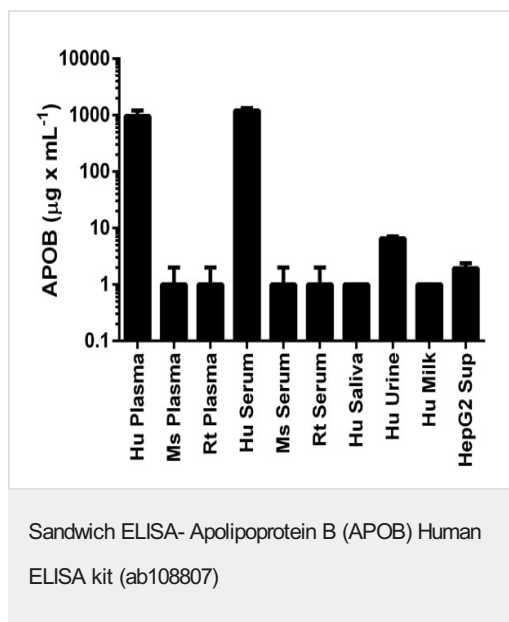
疾病相关 Hypobetalipoproteinemia, familial, 1  
Familial ligand-defective apolipoprotein B-100  
Defects in APOB associated with defects in other genes (polygenic) can contribute to hypocholesterolemia.

序列相似性 Contains 1 vitellogenin domain.

翻译后修饰 Palmitoylated; structural requirement for proper assembly of the hydrophobic core of the lipoprotein particle.

细胞定位 Cytoplasm. Secreted.

## 图片



**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