

### Human Apolipoprotein CI ELISA Kit (APOCI) ab108808

[3 References](#) [1 图像](#)

#### 概述

**产品名称** 人Apolipoprotein CI ELISA试剂盒(APOCI)

**检测方法** Colorimetric

**精确度** 批次内

样品	n	Mean	SD	CV%
Overall				4.8%

批次间

样品	n	Mean	SD	CV%
Overall				9.9%

**样品类型** Cell culture supernatant, Serum, Plasma, Tissue, Cell Lysate

**检测类型** Sandwich (quantitative)

**灵敏度** = 45 ng/ml

**范围** 0.25 µg/ml - 2 µg/ml

**回收率** 98 %

**检测时间** 5h 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 CI (APOCI) Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of apolipoprotein CI concentrations in plasma, serum, cell culture supernatants, cell lysates and tissue samples.

An Apolipoprotein CI specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an Apolipoprotein CI 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 CI 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 M Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human Apolipoprotein C1 Antibody	1 x 120µl
Apolipoprotein C1 Microplate (12 x 8 well strips)	1 unit
Apolipoprotein C1 Standard	2 vials
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml

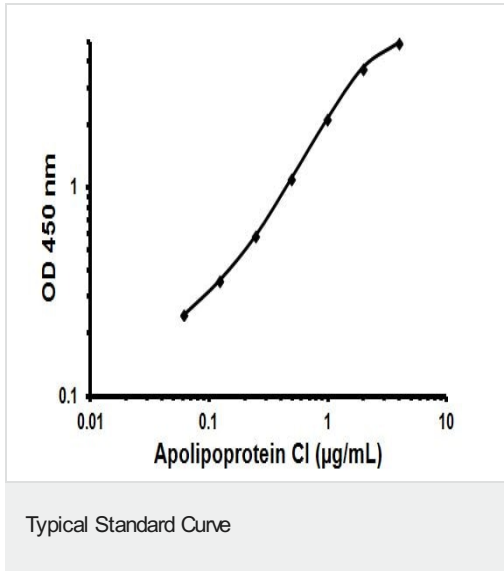
**功能** Appears to modulate the interaction of APOE with beta-migrating VLDL and inhibit binding of beta-VLDL to the LDL receptor-related protein. Binds free fatty acids and reduces their intracellular esterification.

**组织特异性** Synthesized mainly in liver and to a minor degree in intestine. Secreted in plasma.

**序列相似性** Belongs to the apolipoprotein C1 family.

**细胞定位** Secreted.

## 图片



Representative Standard Curve using ab108808

**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