

Human Apolipoprotein CI ELISA Kit (APOCI) ab108808

[3 References](#) [1 Image](#)

Overview

Product name Human Apolipoprotein CI ELISA Kit (APOCI)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				4.8%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				9.9%

Sample type Cell culture supernatant, Serum, Plasma, Tissue, Cell Lysate

Assay type Sandwich (quantitative)

Sensitivity = 45 ng/ml

Range 0.25 µg/ml - 2 µg/ml

Recovery 98 %

Assay time 5h 00m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

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.

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 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

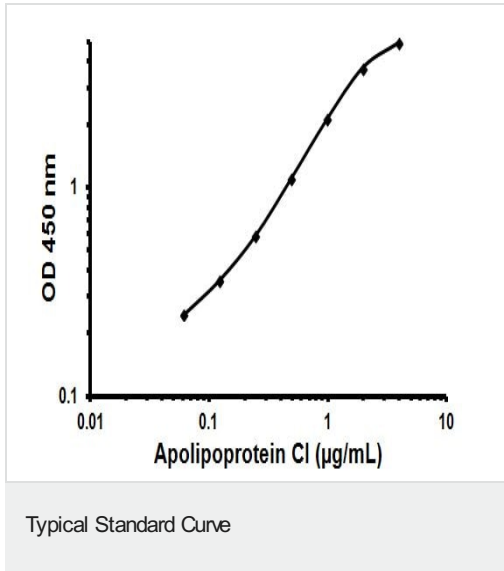
Function Appears to modulate the interaction of APOE with beta-migrating VLDL and inhibit binding of beta-VLDL to the LDL receptor-related protein. Binds free fatty acids and reduces their intracellular esterification.

Tissue specificity Synthesized mainly in liver and to a minor degree in intestine. Secreted in plasma.

Sequence similarities Belongs to the apolipoprotein C1 family.

Cellular localization Secreted.

Images



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.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