



Product datasheet

Human Apolipoprotein CII ELISA Kit ab168549

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

Overview

Product name	Human Apolipoprotein CII ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			6.2%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			10.6%
Sample type	Serum, Plasma			
Assay type	Sandwich (quantitative)			
Sensitivity	= 11 ng/ml			
Range	0.016 µg/ml - 4 µg/ml			
Recovery	98 %			
Assay time	4h 00m			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	Abcam's Apolipoprotein CII Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Apolipoprotein CII concentrations in plasma and serum.			

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

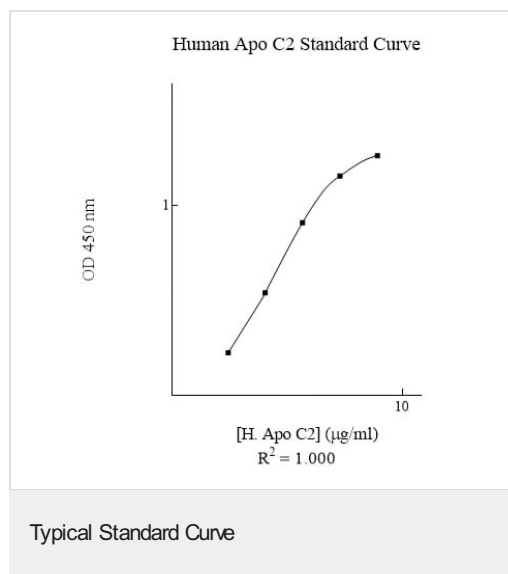
The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

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
40x Biotinylated Human Apolipoprotein CII Antibody	1 x 150µl
Apolipoprotein CII Microplate (12 x 8 well strips)	1 unit
Apolipoprotein CII Standard	2 vials
Chromogen Substrate	1 x 8ml
Sealing Tapes	3 units
Stop Solution	1 x 12ml

Images



Representative Standard Curve using ab168549

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