

Product datasheet

Human Apolipoprotein A-IV Matched Antibody Pair Kit ab220140

Recombinant

2 Images

Overview

Product name	Human Apolipoprotein A-IV Matched Antibody Pair Kit
Detection method	Colorimetric
Assay type	ELISA set
Sensitivity	3.74 pg/ml
Range	250 pg/ml - 16000 pg/ml
Species reactivity	Reacts with: Human
Product overview	<p>Human Apolipoprotein A-IV Matched Antibody Pair Kits include a capture and a biotinylated detector antibody pair, along with a calibrated protein standard, suitable for sandwich ELISA. The Matched Antibody Pair Kit can be used to quantify native and recombinant human Apolipoprotein A-IV.</p> <p>Optimization of the kit reagents to sample type, immunoassay format or instrumentation may be required. Guidelines for use of this kit in a standard 96-well microplate sandwich ELISA using HRP/TMB system of colorimetric detection is described in this assay procedure for the purposes of quantification.</p> <p>Protocol information and tips on the use of the Matched Antibody Pair kits for sandwich ELISA can be found on our website. An accessory pack can be purchased which includes buffer reagents required to perform 10 x 96-well plate sandwich ELISAs (ab210905).</p> <p>For additional information on the performance of the antibody pair used in this kit, please see our equivalent SimpleStep ELISA kit ab214567. Please note that while the antibody pair is the same provided in the corresponding SimpleStep ELISA Kit, due to differences in their formulation, this antibody pair cannot be used with the consumables provided with our SimpleStep ELISA Kits.</p>
Tested applications	Suitable for: ELISA
Platform	Reagents

Properties

Storage instructions

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

Components	10 x 96 tests	5 x 96 tests
Human Apolipoprotein A-IV Capture Antibody	2 x 50µg	1 x 50µg
Human Apolipoprotein A-IV Detector Antibody	2 x 12.5µg	1 x 12.5µg
Human Apolipoprotein A-IV Lyophilized Protein	2 vials	1 vial

Relevance

APOA4 (apolipoprotein A-IV) is a component of HDL and chylomicrons. Its primary site of synthesis is the intestine, in association with lymph chylomicron particles. Although its precise function is not known, APOA4 is a potent activator of lecithin-cholesterol acyltransferase (LCAT) in vitro. In rodents, Apo A-IV inhibits gastric emptying and serves as a satiety factor whose synthesis and secretion are increased by the ingestion of dietary fat. It also possesses anti-inflammatory and antiatherogenic properties

Cellular localization

Secreted

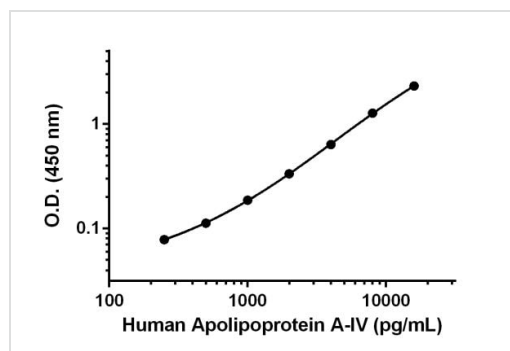
Applications

Our [Abpromise guarantee](#) covers the use of **ab220140** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
ELISA		Use at an assay dependent concentration.

Images



Human Apolipoprotein A-IV standard curve.

Standard calibration curve. Background subtracted values are graphed.

Powered by
recombinant antibodies



Research with confidence
Consistent and reproducible results



Long-term and scalable supply
Recombinant technology



Success from the first experiment
Confirmed specificity



Ethical standards compliant
Animal-free production

Sandwich ELISA - Human Apolipoprotein A-IV Matched Antibody Pair Kit (ab220140)

To learn more about the advantages of recombinant antibodies see [here](#).

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