

Product datasheet

Human Lipoprotein A Matched Antibody Pair Kit ab217612

1 Image

Overview

Product name	Human Lipoprotein A Matched Antibody Pair Kit
Detection method	Colorimetric
Assay type	ELISA set
Range	15.625 ng/ml - 1000 ng/ml
Species reactivity	Reacts with: Human
Product overview	<p>Matched Antibody Pair kit is a titrated unlabeled capture antibody, a titrated biotin-labeled detector and a calibrated protein standard. The Matched Antibody Pair Kit can be used to quantify native and recombinant human Lipoprotein A.</p> <p>Both capture and detector antibodies are rabbit monoclonal antibodies.</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>Additional protocol information and tips on the use of the Matched Antibody Pair kits for sandwich ELISA can be found on our website.</p> <p>For additional information on the performance of the antibody pair used in this kit, please see our equivalent SimpleStep ELISA® (ab212165) which uses the same antibody pair.</p> <p>Buffer information: The capture antibody is glycerol free. The detector antibody contains glycerol.</p> <p>To receive an electronic copy of the Certificate of Analysis, please send an email with "CoA for matched antibody pair kit" in the subject line and the desired product number and lot number in the body of the email.</p>
Tested applications	Suitable for: ELISA
Platform	Reagents

Properties

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

Components	10 x 96 tests	2 x 96 tests
Human Lipoprotein A Capture Antibody	1 x 100µg	1 x 20µg
Human Lipoprotein A Detector Antibody	1 x 25µg	1 x 5µg
Human Lipoprotein A Lyophilized Protein	1 vial	1 vial

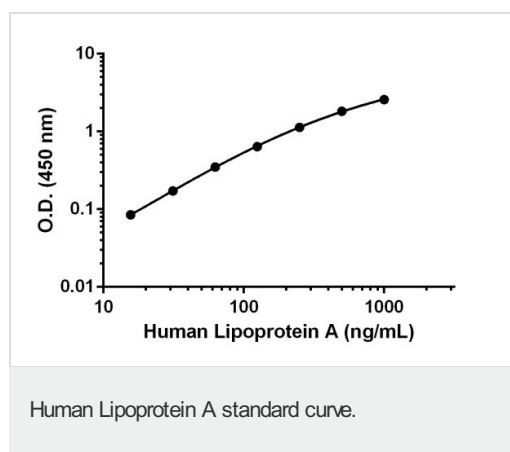
Relevance Lipoprotein(a) (Lp(a)) is a lipoprotein subclass assembled in the blood from low density lipoprotein (LDL) molecules and apolipoprotein-a (apo-a). Lp(a) recruits inflammatory cells through interaction with Mac-1 integrin. High Lp(a) in blood is a risk factor for coronary heart disease, cerebrovascular disease, atherosclerosis, thrombosis, and stroke. Lp(a) concentrations may be affected by disease states, but are only moderately affected by diet, exercise and other environmental factors. Lipid-reducing drugs have no effect on Lp(a) concentration. High Lp(a) predicts risk of early atherosclerosis similar to high LDL, but in advanced atherosclerosis, Lp(a) is a risk factor independent of LDL, indicating a coagulant risk of plaque thrombosis. Apo(a) contains domains that are very similar to plasminogen (PLG). Lp(a) accumulates in the vessel wall and inhibits binding of PLG to the cell surface, reducing plasmin generation which increases clotting. This inhibition also promotes proliferation of smooth muscle cells. These unique features of Lp(a) suggest a role in the generation of clots and atherosclerosis.

Applications

Our [Abpromise guarantee](#) covers the use of **ab217612** 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



Standard calibration curve. Background subtracted values are graphed.

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