

Product datasheet

HDL Uptake Assay Kit (Fluorometric) ab204717

[4 References](#) [2 Images](#)

Overview

Product name	HDL Uptake Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based (quantitative)
Species reactivity	Reacts with: Mammals, Other species
Product overview	HDL Uptake Assay Kit (Fluorometric) (ab204717) contains a fluorescently-labeled HDL that can be measured when taken up by cells. Unlabeled-HDL is included in the kit for assay validation.
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K586 HDL Uptake Assay Kit (Fluorometric). K586-100 is the same size as the 100 test size of ab204717.</p> <p>High-density lipoprotein (HDL) consists of a protein shell, containing apolipoproteins A-I and A-II, and a hydrophobic core consisting of cholesterol and triglycerides surrounded by phospholipids and cholesterol esters. HDL transports cholesterol to the liver or steroidogenic organs such as adrenals, ovary, and testes by both direct and indirect pathways. HDL-Cholesterol is removed by HDL receptors such as scavenger receptor BI (SR-BI), which mediate the selective uptake of cholesterol from HDL. Clinically, high levels of HDL has been associated with cardiovascular health. This is due to the ability of HDL particle to transport cholesterol from lipid-laden macrophages of atherosclerotic arteries to the liver for secretion into the bile by a process called as reverse cholesterol transport.</p>
Platform	Microplate reader

Properties

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

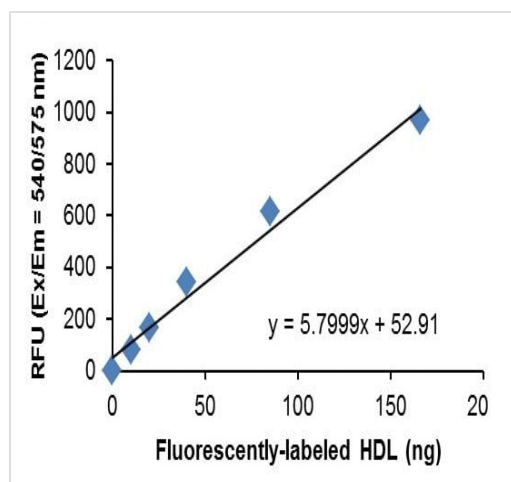
Components	100 tests
Assay Buffer	1 x 100ml
Fluorescently-labeled HDL (5 mg/ml)	1 x 1ml
Unlabeled HDL (2X)	1 x 100µl

Relevance

High-density lipoprotein (HDL) is one of the five major groups of lipoproteins, which, in order of sizes, largest to smallest, are chylomicrons, VLDL, IDL, LDL, and HDL, which enable lipids like cholesterol and triglycerides to be transported within the water-based bloodstream. In healthy individuals, about thirty percent of blood cholesterol is carried by HDL.

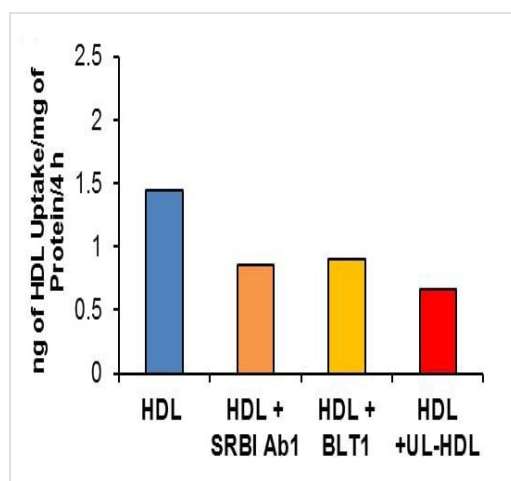
Blood tests typically report HDL-C level, i.e. the amount of cholesterol contained in HDL particles. It is often contrasted with low-density or LDL cholesterol or LDL-C. HDL particles are able to remove cholesterol from within artery atheroma and transport it back to the liver for excretion or re-utilization, which is the main reason why the cholesterol carried within HDL particles (HDL-C) is sometimes called "good cholesterol" (despite the fact that it is exactly the same as the cholesterol in LDL particles).

Images



Fluorescently-labeled HDL Standard curve.

HDL Uptake Assay Kit (Fluorometric) (ab204717)



Fluorescently-labeled HDL uptake in HepG2 (hepatocytes) cells. Cells were starved for 4 hrs and incubated with Fluorescently-labeled HDL (10 μ g) for 4 hrs. Fluorescently-labeled HDL (HDL) uptake was inhibited by SR-BI (HDL receptor) antibody (HDL + SRB1 Ab1), BLT1 (SR-BI inhibitor; HDL + BLT1) and outcompeted by Unlabeled-HDL (HDL+ UL-HDL [20 μ g]). Background was subtracted from the data (Wash off was subtracted as background).

HDL Uptake Assay Kit (Fluorometric) (ab204717)

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