

Product datasheet

Cholesterol Assay Kit (Cell-Based) ab133116

★★★★★ 1 Abreviews 19 References 6 Images

Overview

Product name	Cholesterol Assay Kit (Cell-Based)
Detection method	Fluorescent
Sample type	Adherent cells
Assay type	Cell-based
Assay time	1h 15m
Species reactivity	Reacts with: Mammals, Other species
Product overview	Cholesterol Assay Kit (Cell-Based) (ab133116) includes filipin III, fixative, and wash buffer in a ready to use format. It provides a simple fluorometric method to study mechanisms and biological factors that regulate cholesterol metabolism or movement within cells. A cholesterol trafficking inhibitor, U-18666A, is included as a positive control.

Cholesterol assay protocol summary:

- remove medium from cells
- fix with fixative solution for 10 min
- wash with wash buffer 3 times for 5 min
- add Filipin III solution and incubate for 30-60 min
- wash twice for 5 min
- analyze with fluorescent microscope

Notes Cholesterol is both an important structural component of cell membranes and an early intermediate in hormone and bile acid biosynthesis. Cholesterol is not uniformly distributed among cellular membranes / organelle membranes, but rather there are structurally and kinetically distinct cholesterol rich and poor domains.

Filipin III is the predominant isomer of filipin, the collective name given to four isomeric polyene macrolides isolated from cultures of *S. filipinensis*. Filipin is widely used as a probe for sterol location in biological membranes.

Interaction with cholesterol alters the filipin absorption and fluorescence spectra allowing visualization with a fluorescence microscope capable of excitation at 340-380 nm and emission at 385-470 nm. Filipin's ease of use makes it a convenient tool for the histochemical identification of unesterified cholesterol both *in vitro* and *in vivo*.

Other cholesterol assay kits include:

- [HDL and LDL/LDL Cholesterol assay kit ab65390](#)
- [Cholesterol/Cholesterol Ester assay kit ab65359](#)
- [Cholesterol Efflux assay kit ab196985](#)

- [Cholesterol Uptake assay kit ab236212](#)

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

Platform Fluorescence microscope

Properties

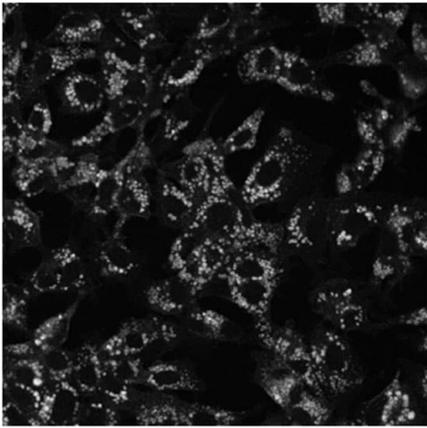
Storage instructions Please refer to protocols.

Components	192 tests
Cell-Based Assay Fixative	2 vials
Cell-Based Assay U-18666A	1 vial
Cholesterol Detection Assay Buffer	1 vial
Cholesterol Detection Filipin III	1 vial
Cholesterol Detection Wash Buffer	1 vial

Relevance

Cholesterol is a waxy substance that occurs naturally in all parts of the body and that the body needs to function normally. It is present in cell walls or membranes everywhere in the body, including the brain, nerves, muscle, skin, liver, intestines, and heart. The body uses cholesterol to produce many hormones, vitamin D, and the bile acids that help to digest fat. It takes only a small amount of cholesterol in the blood to meet these needs. However, if you have too much cholesterol in your bloodstream, it can lead to atherosclerosis, a condition in which fat and cholesterol are deposited in the walls of the arteries in many parts of the body, including the coronary arteries feeding the heart. In time, narrowing of the coronary arteries by atherosclerosis can produce the signs and symptoms of heart disease, including angina and heart attack.

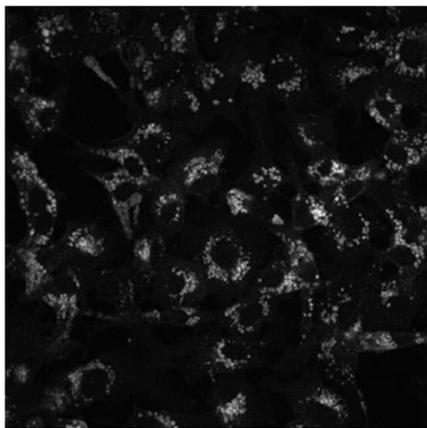
Images



Cholesterol Assay Kit (Cell-Based) (ab133116)

HepG2 cells in response to 2 μM U-18666A.

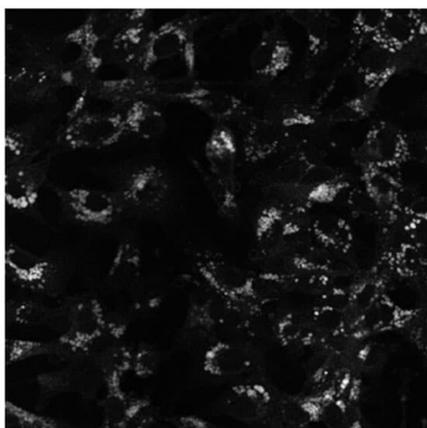
HepG2 cells were seeded in a 96-well plate at a density of 2.4×10^4 cells/well and cultured overnight. The next day, cells were treated with 2 μM U-18666A for 48 hours. U-18666A treatment for 48 hours induces intracellular accumulation.



Cholesterol Assay Kit (Cell-Based) (ab133116)

HepG2 cells in response to 1 μM U-18666A.

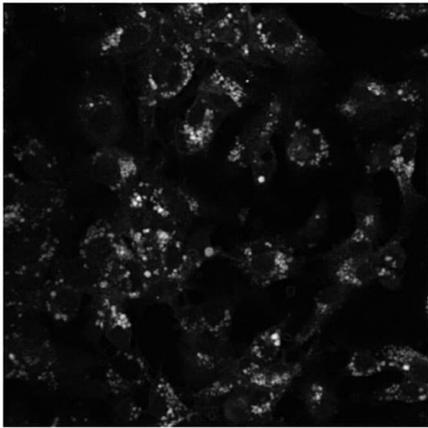
HepG2 cells were seeded in a 96-well plate at a density of 2.4×10^4 cells/well and cultured overnight. The next day, cells were treated with 1 μM U-18666A for 48 hours. U-18666A treatment for 48 hours induces intracellular accumulation.



Cholesterol Assay Kit (Cell-Based) (ab133116)

HepG2 cells in response to 0.5 μM U-18666A.

HepG2 cells were seeded in a 96-well plate at a density of 2.4×10^4 cells/well and cultured overnight. The next day, cells were treated with 0.5 μM U-18666A for 48 hours. U-18666A treatment for 48 hours induces intracellular accumulation.

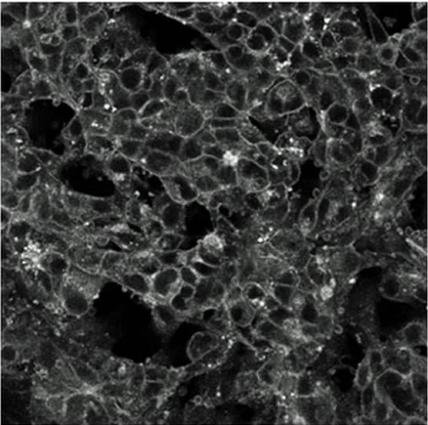


Copyright (c) 2015 Abcam Plc

Cholesterol Assay Kit (Cell-Based) (ab133116)

HepG2 cells in response to 0.25 μM U-18666A.

HepG2 cells were seeded in a 96-well plate at a density of 2.4×10^4 cells/well and cultured overnight. The next day, cells were treated with 0.25 μM U-18666A for 48 hours. U-18666A treatment for 48 hours induces intracellular accumulation.

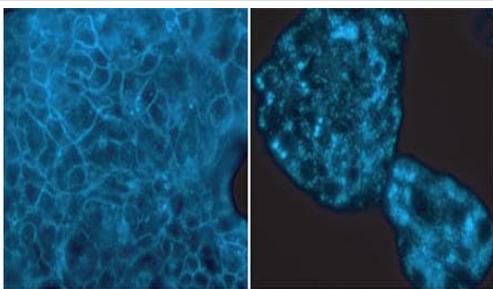


Copyright (c) 2015 Abcam Plc

Cholesterol Assay Kit (Cell-Based) (ab133116)

Accumulation of cholesterol inside HepG2 cells - Control.

HepG2 cells were seeded in a 96-well plate at a density of 2.4×10^4 cells/well and cultured overnight. The next day, cells were treated with DMSO (vehicle) for 48 hours. Cells demonstrate that majority of cholesterol is localized on the plasma membrane.



Cholesterol Assay Kit (Cell-Based) (ab133116)

Accumulation of cholesterol inside HepG2 cells in response to 1.25 μM U-18666A.

HepG2 cells were seeded in a 96-well plate at a density of 3×10^4 cells/well and cultured overnight. The next day, cells were treated with DMSO (vehicle) or 1.25 μM U-18666A for 48 hours. Panel A: Cells treated with DMSO alone demonstrate that majority of cholesterol is localized on the plasma membrane. Panel B: U-18666A treatment for 48 hours induces intracellular accumulation.

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