

LDL Uptake Assay Kit (Cell-Based) ab133127

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

Overview

Product name	LDL Uptake Assay Kit (Cell-Based)
Detection method	Fluorescent
Sample type	Adherent cells
Assay type	Cell-based
Assay time	6h 30m
Species reactivity	Reacts with: Mammals, Other species
Product overview	<p>LDL Uptake Assay Kit (Cell-Based) (ab133127) provides a convenient tool for studying LDL uptake and regulation at the cellular level. This kit employs Human LDL conjugated to DyLight™ 550 as a fluorescent probe for detection of LDL uptake into cultured cells. An LDL receptor-specific polyclonal antibody and a DyLight™ 488-conjugated secondary antibody are included for identifying the distribution of LDL receptors.</p> <p>LDL uptake assay protocol summary:</p> <ul style="list-style-type: none"> - remove culture medium from experimentally treated cells - add LDL-Dylight 550 solution and incubate for 3-4 hrs - replace solution with culture medium - analyze LDL uptake with fluorescent microscope - wash cells and fix for 10 min - wash 3 times for 5 min - add assay blocking solution and incubate for 30 min - add anti-LDL receptor antibody and incubate for 1 hr - wash 3 times for 5 min - add Dylight 488 secondary antibody and incubate for 1 hr - wash 3 times for 5 min - analyze staining with fluorescence microscope
Notes	<p>LDL is the major carrier of cholesterol in the blood, accounting for more than 60% of total plasma cholesterol. LDL is taken up by hepatic and extra-hepatic tissues through receptor mediated endocytosis triggered by apoB-100-LDL receptor interaction. The internalized LDL particle is transported to lysosomes where it is degraded to free cholesterol and amino acids.</p>
Platform	Fluorescence microscope

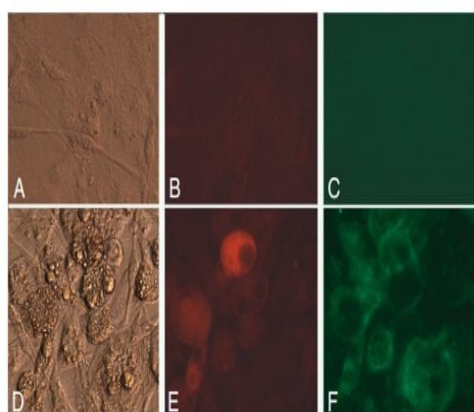
Properties

Storage instructions

Please refer to protocols.

Components	1 kit
Cell-Based Assay Blocking Solution	1 vial
Cell-Based Assay Fixative	1 vial
DyLight 488-Conjugated Goat Anti-Rabbit Secondary Antibody	1 unit
LDL-DyLight 550	1 vial
Rabbit Anti-LDL Receptor Primary Antibody	1 vial

Images



Functional Studies - LDL Uptake Assay Kit (Cell-Based) (ab133127)

LDL uptake in pre-adipocytes and adipocytes.

Panel A: Light image of undifferentiated pre-adipocytes.

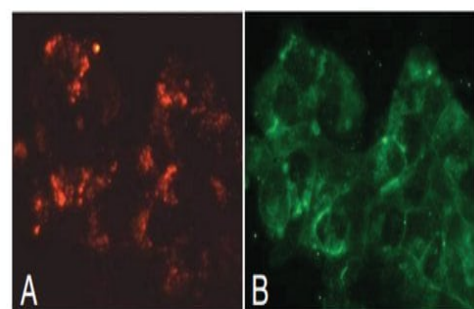
Panel B: Undifferentiated pre-adipocytes do not show any LDL uptake.

Panel C: Undifferentiated pre-adipocytes show little LDL receptor expression.

Panel D: Light image of differentiated adipocytes.

Panel E: Corresponding cells to those in Panel D showing uptake of LDL (red).

Panel F: Corresponding cells to those in Panel D showing expression of LDL receptor (green).



LDL Uptake Assay Kit (Cell-Based) (ab133127)

LDL uptake in HepG2 cells.

HepG2 cells were cultured at a density of 3×10^4 cells/well in a 96 well plate for two days then treated with 32.5 μ M EGCG overnight. LDL-Dylight™ 549 (10 μ g/ml) was added and the cells were incubated for an additional four hours. Cells were fixed and stained for LDL receptor using a Rabbit anti-LDL receptor primary antibody and DyLight™ 488-conjugated secondary antibody.

Panel A: LDL-Dylight™ 549 taken into cells appears in red.

Panel B: LDL receptors (in green) show a distribution pattern that matches the cells in the Panel A containing LDL-Dylight™ 549.

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