

Product datasheet

Cholesterol Uptake Assay Kit ab236212

[2 References](#) [3 Images](#)

Overview

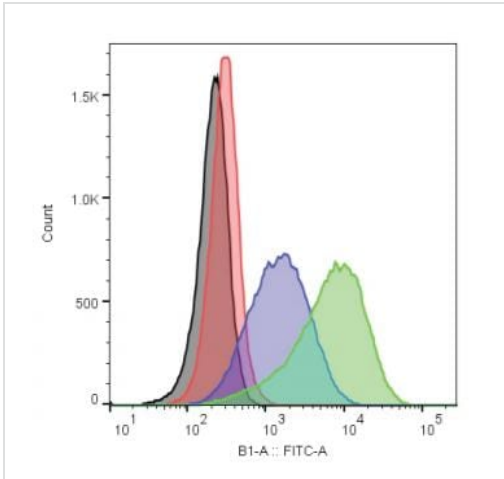
Product name	Cholesterol Uptake Assay Kit
Detection method	Fluorescent
Sample type	Adherent cells, Suspension cells
Product overview	Cholesterol Uptake Assay Kit (ab236212) provides a convenient tool for studying cellular cholesterol trafficking. The kit employs NBD Cholesterol, a fluorescently-tagged cholesterol, as a probe for the detection of cholesterol taken up by cultured cells. U-18666A, which increases cholesterol uptake by inhibiting trafficking of synthesized cholesterol, is included as a positive control. The kit provides enough NBD Cholesterol to test 250 samples in a 96-well format.
Notes	Other cholesterol assay kits include: <ul style="list-style-type: none"> - HDL and LDL/LDL Cholesterol assay kit ab65390 - Cell-based Cholesterol assay kit ab133116 - Cholesterol/Cholesterol Ester assay kit ab65359 - Cholesterol Efflux assay kit ab196985
Platform	Microplate reader, Fluor. microscope, Flow cyt.

Properties

Storage instructions Please refer to protocols.

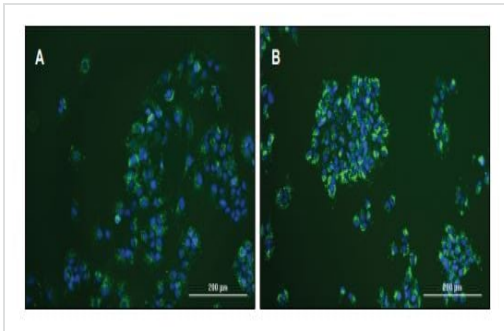
Components	1 kit
Cell-Based Assay Buffer Tablet	1 tablet
Cell-Based Assay NBD Cholesterol	1 x 500µl
Cell-Based Assay U-18666A	1 x 100µl

Images

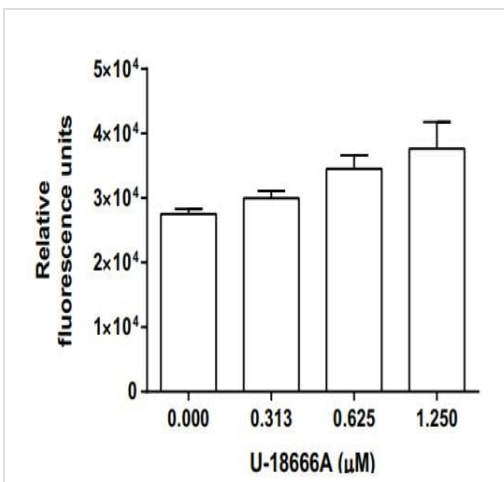


U-18666A increases cholesterol uptake in Jurkat cells as measured by flow cytometry

Jurkat cells were seeded at a density of 5×10^5 cells/mL and incubated overnight in serum-free RPMI with U-18666A or vehicle and 20 $\mu\text{g/mL}$ NBD Cholesterol in a cell culture incubator at 37°C . The next day, cells were transferred to a v-bottom plate for washing and reading on a flow cytometer. Cholesterol uptake was evaluated in the live cell gate using FlowJo analysis software. U-18666A at both 2.5 μM (green) and 1.25 μM (blue) showed a significant ($p < 0.05$, t-test) shift in mean fluorescence as compared to the vehicle control (black) and the untreated (red) cells.



Blocking intracellular cholesterol transport with U-18666A increases NBD cholesterol uptake



U-18666A causes a dose-dependent increase in NBD Cholesterol uptake in Caco-2 cells, as measured on a fluorescent plate reader

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