

# Cholesterol Uptake Assay Kit ab236212

**3 References**   [画像数 3](#)

### 製品の概要

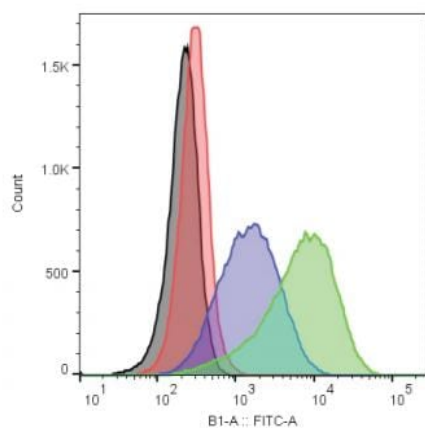
製品名	Cholesterol Uptake Assay Kit
検出方法	Fluorescent
サンプルの種類	Adherent cells, Suspension cells
製品の概要	Cholesterol Uptake Assay Kit (ab236212) provides a convenient tool for studying cellular cholesterol trafficking. The kit employs NBD Cholesterol, a fluorescently-tagged cholesterol, as a probe for the detection 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.
特記事項	Other cholesterol assay kits include: - <b><u>HDL and LDL/VLDL Cholesterol assay kit ab65390</u></b> - <b><u>Cell-based Cholesterol assay kit ab133116</u></b> - <b><u>Cholesterol/Cholesterol Ester assay kit ab65359</u></b> - <b><u>Cholesterol Efflux assay kit ab196985</u></b>
試験プラットフォーム	Microplate reader, Fluor. microscope, Flow cyt.

### 製品の特性

**保存方法** Please refer to protocols.

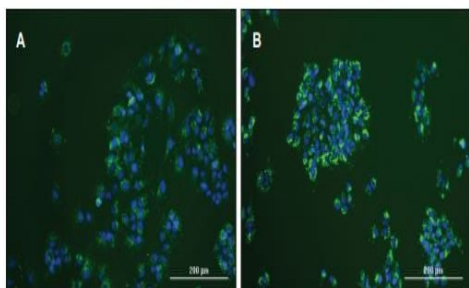
内容	1 kit	1 kit
Cell-Based Assay Buffer Tablet	1 tablet	1 tablet
Cell-Based Assay NBD Cholesterol	1 x 500µl	1 x 500µl
Cell-Based Assay U-18666A	1 x 100µl	1 x 100µl

### 画像

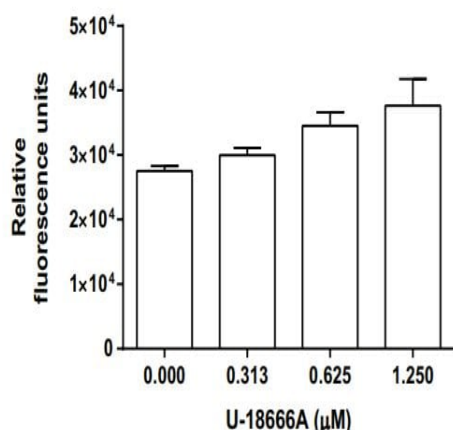


U-18666A increases cholesterol uptake in Jurkat cells as measured by flow cytometry

Jurkat cells were seeded at a density of  $5 \times 10^5$  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  $37^\circ\text{C}$ . 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  $\mu\text{M}$  (green) and 1.25  $\mu\text{M}$  (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.co.jp/abpromise> or contact our technical team.

### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors