

Product datasheet

Cholesterol Efflux Assay Kit (Cell-based) ab196985

1 References [画像数 1](#)

製品の概要

<b>製品名</b>	Cholesterol Efflux Assay Kit (Cell-based)
<b>検出方法</b>	Fluorescent
<b>サンプルの種類</b>	Adherent cells, Purified protein
<b>アッセイタイプ</b>	Cell-based (quantitative)
<b>種交差性</b>	<b>交差種:</b> Mammals, Other species
<b>製品の概要</b>	Cholesterol Efflux Assay Kit ab196985 is a high-throughput, cell-based screening assay for measuring cholesterol efflux. The cholesterol efflux assay protocol uses fluorescently-labeled cholesterol (Ex/Em 482/515 nm). It provides a safe, sensitive and reproducible method for measuring cholesterol efflux.

This cholesterol efflux assay can be used to:  
 - screen small molecules for their effect on cholesterol efflux as a part of drug discovery program.

Cholesterol efflux assay protocol summary:  
 - label cells with labeling and equilibration mix overnight  
 - wash cells, treat as required and incubate  
 - transfer cell supernatant to microplate, and separately solubilize cells with cell lysis buffer  
 - analyze supernatant and cell lysates with microplate reader

**特記事項**  
 Cholesterol efflux from the peripheral tissues and cells in atherosclerotic plaque is an initial and critical step in Reverse Cholesterol Transport (RCT). RCT is the process by which extra hepatic cells, including macrophage-derived foam cells in arterial atherosclerotic plaque, transport excessive cholesterol back to the liver for bile acid synthesis and excretion, thus lowering the peripheral lipid burden.

**試験プラットフォーム**                      Microplate reader

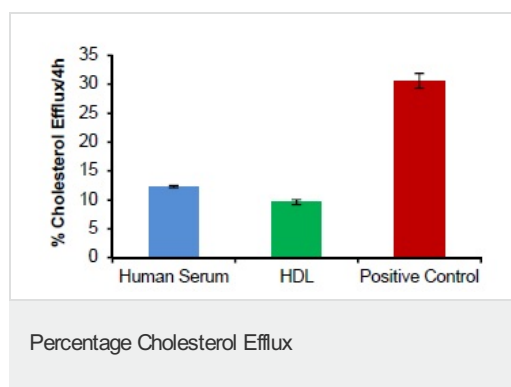
製品の特性

**保存方法**                                      Store at -20°C. Please refer to protocols.

内容	100 tests
Cell Lysis Buffer	1 x 20ml

内容	100 tests
Equilibration Buffer	1 x 5ml
Labeling Reagent	1 x 5ml
Positive Control	1 x 1ml
Reagent A	1 x 10 $\mu$ l
Reagent B	1 x 10 $\mu$ l
Serum Treatment Reagent	1 x 1ml

## 画像



J774.1 cells were labeled with the Labeling Media and treated with various cholesterol acceptors like Human Serum, HDL (50  $\mu$ g) or Positive control known to cause cholesterol efflux. Cholesterol efflux is expressed as % efflux elicited by cells in 4 hours.

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

## 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