

Cholesterol Efflux Assay Kit (Cell-based) ab196985

30 References **1 图像**

概述

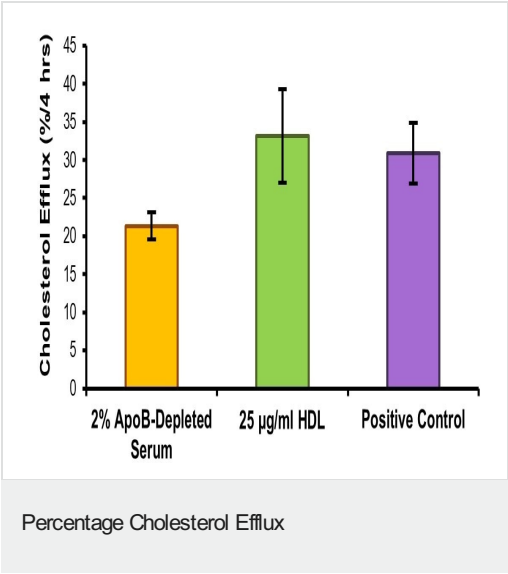
产品名称	Cholesterol Efflux Assay试剂盒(Cell-based)
检测方法	Fluorescent
样品类型	Adherent cells, Purified protein
检测类型	Cell-based (quantitative)
种属反应性	与反应: Mammals, Other species
产品概述	<p>Cholesterol Efflux Assay Kit ab196985 is a high-throughput, cell-based screening assay for measuring cholesterol efflux. The cholesterol efflux assay protocol uses fluorescently-labeled cholesterol (Ex/Em 482/515 nm). It provides a safe, sensitive and reproducible method for measuring cholesterol efflux.</p> <p>This cholesterol efflux assay can be used to:</p> <ul style="list-style-type: none"> - screen small molecules for their effect on cholesterol efflux as a part of drug discovery program. <p>Cholesterol efflux assay protocol summary:</p> <ul style="list-style-type: none"> - 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
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K582 Cholesterol Efflux Fluorometric Assay Kit (cell-based). K582-100 is the same size as the 100 test size of ab196985.</p> <p>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.</p> <p>Other cholesterol assay kits include:</p> <ul style="list-style-type: none"> - <u>HDL and LDL/VLDL Cholesterol assay kit ab65390</u> - <u>Cell-based Cholesterol assay kit ab133116</u> - <u>Cholesterol/Cholesterol Ester assay kit ab65359</u> - <u>Cholesterol Uptake assay kit ab236212</u>
平台	Microplate reader

性能

存放说明 Store at -20°C. Please refer to protocols.

组件	100 tests
Cholesterol Uptake Positive Control	1 x 1ml
Equilibration Buffer	1 x 5ml
Labeling Reagent	1 x 5ml
Lysis Buffer I	1 x 25ml
Reagent A	1 x 10µl
Reagent B	1 x 50µl
Serum Treatment Reagent	1 x 1ml

图片



Percentage (%) Labeled Cholesterol Efflux. J774.1 cells were labeled with the fluorescent cholesterol Labeling Media and treated with various cholesterol acceptors such as 2% LDL/VLDL-depleted Human Serum (2 µl/well), purified human HDL (25 µg/ml, 2.5 µg/well) or Positive Control reagent. Cholesterol efflux is expressed as percentage of fluorescently-labeled cholesterol transferred from labeled adherent macrophages to designated cholesterol acceptor present in supernatant over a 4 hour time period, with no treatment control (efflux to medium containing no acceptor) percentage subtracted (mean ± SD of N = 5-6 wells).

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Promise 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.cn/abpromise> or contact our technical team.

Terms and conditions

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