

Free Fatty Acid Uptake Assay Kit (Fluorometric) ab176768

★☆☆☆☆ [1 Abreviews](#) [21 References](#) [2 图像](#)

概述

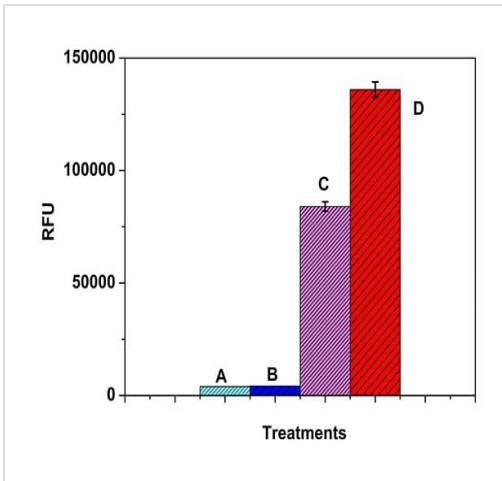
产品名称	Free Fatty Acid Uptake Assay试剂盒(Fluorometric)
检测方法	Fluorescent
样品类型	Adherent cells
检测类型	Cell-based (quantitative)
检测时间	1h 00m
种属反应性	与反应: Mammals, Other species
产品概述	Free Fatty Acid Uptake Assay Kit (Fluorometric) (ab176768) provides a simple and sensitive method for the measurement of fatty acid uptake in cells containing fatty acid transporters. The kit uses a proprietary dodecanoic acid fluorescent fatty acid substrate. This fatty acid uptake assay kit can be performed on any fluorescence microplate reader with a bottom-read mode at Ex/Em = 485/515 nm or FITC channel. The assay can be performed in 96-well or 384-well microtiter plates in a simple mix-and-read procedure, and easily adapted for high throughput screening applications.
说明	Fatty acid uptake is an important therapeutic target for the treatment of many human diseases such as obesity, type 2 diabetes and hepatic steatosis.
平台	Microplate reader

性能

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

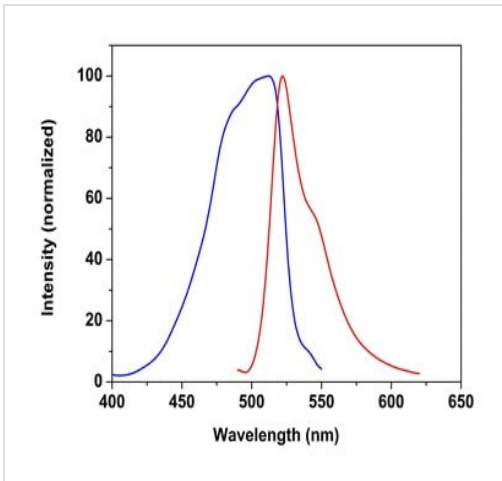
组件	100 tests
Assay Buffer	1 x 10ml
DMSO	1 x 100µl
TF2-C12 Fatty Acid	1 vial

图片



Comparison of fatty acid uptake by 3T3-L1 adipocytes and fibroblast.

Cells were plated at 50,000 cells/100 μ L/well in a 96 well black wall/clear bottom poly-D lysine plate for 5 hours, and then serum deprived for 1 hour. Cells were treated without (control) or with insulin (150 nM), and incubated at 37 $^{\circ}$ C, 5% CO₂ incubator for 30 min. At the end of the incubation time, 100 μ L of fatty acid mixture was added into the well, and incubated for another 60 min, the fluorescence signal was measured with a plate reader using bottom read mode. A – fibroblasts (Control); B – fibroblasts (Insulin); C – adipocytes (Control); D– adipocytes (Insulin).



Excitation and Emission Spectra of Free Fatty Acid Uptake Assay Kit (Fluorometric) (ab176768)

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.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