

**ab176768**

**Free Fatty Acid Uptake  
Assay Kit (Fluorometric)**

**Instructions for Use**

For measurement of fatty acid uptake in cells containing fatty acid transporters.

This product is for research use only and is not intended for diagnostic use.

Version: 2b Last Updated: 8 January 2021

# Table of Contents

---

Table of Contents	2
1. Overview	3
2. Protocol Summary	4
3. Kit Components	5
4. Storage and Stability	5
5. Materials Required, Not Supplied	6
6. Assay Protocol	7
7. Data Analysis	10
8. Troubleshooting	13

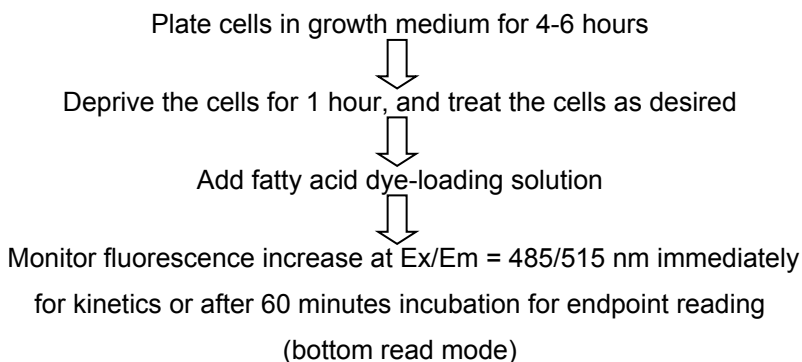
# 1. Overview

---

Fatty acid uptake is an important therapeutic target for the treatment of many human diseases such as obesity, type 2 diabetes and hepatic steatosis. Abcam's 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.

## 2. Protocol Summary

---



## 3. Kit Components

---

Item	Quantity	Storage upon arrival	Storage after use/ reconstitution
TF2-C12 Fatty Acid	1 vial	-20°C	-20°C
Assay Buffer	10 mL	-20°C	-20°C
DMSO	100 µL	-20°C	-20°C

## **4. Storage and Stability**

---

Upon arrival, store the kit at -20°C and protected from light. Please read the entire protocol before performing the assay. Avoid repeated freeze/thaw cycles.

Warm all buffers to room temperature before use. Briefly centrifuge all small vials prior to opening.

## **5. Materials Required, Not Supplied**

---

- 96 or 384-well black plate with clear flat bottoms
- 75cm flask
- Serum free medium
- Fluorescence microplate reader

## 6. Assay Protocol

---

### 1. Reagent Preparation

Thaw all the kit components at room temperature before use.

#### a) **TF2-C12 Fatty Acid stock solution:**

Add 20  $\mu$ L DMSO into the vial of the TF2-C12 Fatty Acid, and mix them well.

*Note: 20  $\mu$ L of the fluorescent fatty acid substrate stock solution is enough for one plate. The unused fluorescent fatty acid substrate stock solution can be aliquoted and stored at  $< -20$  °C for up to two months if the tubes are sealed tightly and kept from light. Avoid repeated freeze-thaw cycles.*

#### b) **Fatty acid dye-loading solution:**

Add 20  $\mu$ L of the TF2-C12 Fatty Acid stock solution (from Step 1a) to 10 mL of Assay Buffer, and mix them well.

*Note: 10 mL of fatty acid dye-loading solution is enough for one plate, prepare fresh for each experiment.*

### 2. Prepare cells

Prepare cells as desired. The following protocols are guidelines to prepare 3T3-L1 adipocytes.

#### a) Prepare differentiated 3T3-L1 adipocytes: 3T3-L1 cells were grown 2 days in a 75 cm flask post-confluence in

DMEM/FBS, and then for 2 days more in DMEM/FBS supplemented with 0.83 $\mu$ M insulin, 0.25 $\mu$ M dexamethasone, and 0.25mM isobutylmethylxanthine.

The medium is changed to maintain the insulin concentration with dexamethasone and IBMX absent for another 2 days. The medium was then changed to DMEM/FBS alone for another 3-5 days.

Differentiated cells (at least 95% of which showed an adipocyte phenotype by accumulation of lipid droplets) were used on day 8 to 12 after induction of differentiation.

- b)** Plate 3T3-L1 adipocytes in growth medium at  $5 \times 10^4$  –  $8 \times 10^4$  cells/well/100 $\mu$ L/ 96-well or 12,500-20,000 cells/well/25 $\mu$ L/ 384-well black wall or clear bottom cell culture Poly-D lysine plate for 4-6 hours before experiment. Centrifuge the plate at 800 rpm for 2 minutes with brake off.

*Note 1: It is recommended to plate 3 wells with growth medium only (without cells) as blank wells for data normalization;*

*Note 2: We find that adipocytes plated at the same day (4-6 hours, and then serum deprived for 1 hour) give better results than that plated for overnight.*

- c)** Remove the cell plate from the incubator, aspirate the medium from the wells, and deprive the cells with 90 $\mu$ L/well/ 96 well-plate or 20  $\mu$ L/well/384 well-plate serum free medium. Incubate the cells at 37 °C, 5% CO<sub>2</sub> incubator for 1 hr.

- d)** Treat the cells by adding 10 $\mu$ L/well/96-well plate (5 $\mu$ L/well/384-well plate) of the test compounds or 1X Hanks and 20 mM Hepes buffer (1X HBSS, pH 7.4) or buffer of your choice as the compound diluent. For blank wells, add the compound diluents. Incubate the cells at 37°C, 5% CO<sub>2</sub> incubator for a desired period of time (30 minutes for 3T3-L1 cells treated with Insulin).

### **3. Run Fatty Acid Uptake assay**

- a)** Remove compound-treated cell plates from the incubator (from Step 2d)' add 100 $\mu$ L/well/96-well plate or 25 $\mu$ L/well/384-well plate (including blank wells) of the fatty acid dye-loading solution (from Step 1b).
- b)** Measure the fluorescence signal with a fluorescence microplate reader at Ex/Em = 485/515 nm (cut off at 495 nm) using a bottom read mode.

For kinetic reading: Read the fluorescence intensity immediately at 20 seconds interval for 30-60 minutes.

For endpoint reading: Read the fluorescence intensity at the end of the 30-60 minutes incubation.

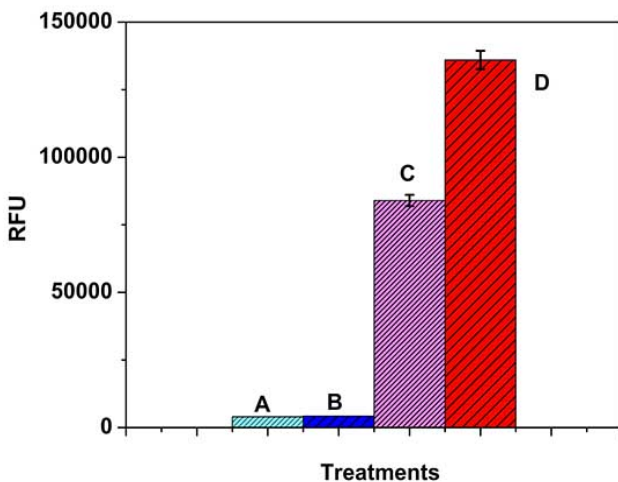


## 7. Data Analysis

---

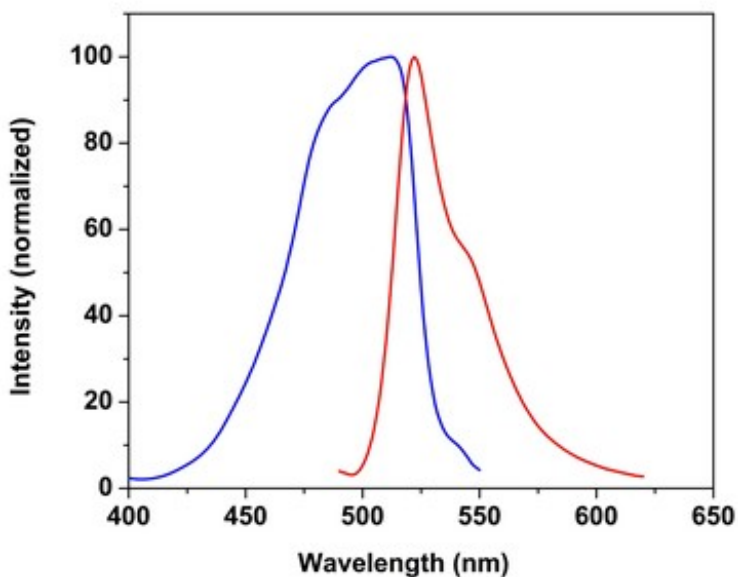
### Calculations:

The fluorescence in blank wells with the growth medium is subtracted from the values for the wells with cells. The background fluorescence of the blank wells may vary depending on the sources of the growth media or the microtiter plates.



**Figure 1.** Comparison of fatty acid uptake by 3T3-L1 adipocytes and fibroblast. Cells were plated at 50,000 cells/100  $\mu$ 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 °C, 5% CO<sub>2</sub> incubator for 30 min. At the end of the incubation time, 100 μL 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).



**Figure 2.** Excitation and emission spectra of Free Fatty Acid Uptake Assay Kit (Fluorometric) (ab176768).

## 8. Troubleshooting

<b>Problem</b>	<b>Reason</b>	<b>Solution</b>
Assay not working	Assay buffer at wrong temperature	Assay buffer must not be chilled - needs to be at RT
	Protocol step missed	Re-read and follow the protocol exactly
	Plate read at incorrect wavelength	Ensure you are using appropriate reader and filter settings (refer to datasheet)
	Unsuitable microtiter plate for assay	Fluorescence: Black plates (clear bottoms); Luminescence: White plates; Colorimetry: Clear plates. If critical, datasheet will indicate whether to use flat- or U-shaped wells
Unexpected results	Measured at wrong wavelength	Use appropriate reader and filter settings described in datasheet
	Samples contain impeding substances	Troubleshoot and also consider deproteinizing samples
	Unsuitable sample type	Use recommended samples types as listed on the datasheet
	Sample readings are outside linear range	Concentrate/ dilute samples to be in linear range

<b>Problem</b>	<b>Reason</b>	<b>Solution</b>
Samples with inconsistent readings	Unsuitable sample type	Refer to datasheet for details about incompatible samples
	Samples prepared in the wrong buffer	Use the assay buffer provided (or refer to datasheet for instructions)
	Samples not deproteinized (if indicated on datasheet)	Use the <b>10kDa spin column (ab93349)</b>
	Cell/ tissue samples not sufficiently homogenized	Increase sonication time/ number of strokes with the Dounce homogenizer
	Too many freeze-thaw cycles	Aliquot samples to reduce the number of freeze-thaw cycles
	Samples contain impeding substances	Troubleshoot and also consider deproteinizing samples
	Samples are too old or incorrectly stored	Use freshly made samples and store at recommended temperature until use
Lower/ Higher readings in samples and standards	Not fully thawed kit components	Wait for components to thaw completely and gently mix prior use
	Out-of-date kit or incorrectly stored reagents	Always check expiry date and store kit components as recommended on the datasheet
	Reagents sitting for extended periods on ice	Try to prepare a fresh reaction mix prior to each use
	Incorrect incubation time/ temperature	Refer to datasheet for recommended incubation time and/ or temperature
	Incorrect amounts used	Check pipette is calibrated correctly (always use smallest volume pipette that can pipette entire volume)

<b>Problem</b>	<b>Reason</b>	<b>Solution</b>
Standard curve is not linear	Not fully thawed kit components	Wait for components to thaw completely and gently mix prior use
	Pipetting errors when setting up the standard curve	Try not to pipette too small volumes
	Incorrect pipetting when preparing the reaction mix	Always prepare a master mix
	Air bubbles in wells	Air bubbles will interfere with readings; try to avoid producing air bubbles and always remove bubbles prior to reading plates
	Concentration of standard stock incorrect	Recheck datasheet for recommended concentrations of standard stocks
	Errors in standard curve calculations	Refer to datasheet and re-check the calculations
	Use of other reagents than those provided with the kit	Use fresh components from the same kit

## Notes



**For all technical and commercial enquires please go to:**

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)