

Screen Quest™ Fluorimetric Fatty Acid Uptake Assay Kit

Catalog number: 36385
Unit size: 100 Tests

Component	Storage	Amount
Component A: TF2-C12 Fatty Acid	Freeze (<-15 °C), Minimize light exposure	1 vial, lyophilized
Component B: Assay Buffer	Freeze (<-15 °C), Minimize light exposure	10 mL
Component C: DMSO	Freeze (<-15 °C)	100 µL

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. The Screen Quest™ Fluorimetric Fatty Acid Uptake Assay Kit 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.

AT A GLANCE

Protocol summary

1. Plate cells in growth medium for 4-6 hours
2. Transfer the cells into serum free medium for 1 hour and treat cells as desired
3. Add 100 µL/well of the Fatty Acid dye-loading solution
4. Monitor fluorescence increase at Ex/Em = 485/515 nm immediately for kinetics or after 60 minutes incubation for endpoint reading (bottom read mode)

Important Thaw all the kit components at room temperature before starting the experiment.

KEY PARAMETERS

Instrument:	Fluorescence microplate reader
Excitation:	485 nm
Emission:	515 nm
Cutoff:	495 nm
Recommended plate:	Solid black

PREPARATION OF STOCK SOLUTIONS

Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20 °C after preparation. Avoid repeated freeze-thaw cycles.

TF2-C12 Fatty Acid stock solution:

Add 20 µL of DMSO (Component C) to the vial of TF2-C12 Fatty Acid (Component A) and mix them well.

Note 20 µL 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

PREPARATION OF WORKING SOLUTION

Fatty Acid dye-loading solution:

Add 20 µL of the TF2-C12 Fatty Acid stock solution to 10 mL of Assay Buffer (Component B) and mix them well.

Note 10 mL of Fatty Acid dye-loading solution is enough for one plate; prepare fresh for each plate and experiment.

SAMPLE EXPERIMENTAL PROTOCOL

1. Treat cells with test compounds as desired.
2. Remove compound-treated cell plates from the incubator, add 100 µL/well (96-well plate) or 25 µL/well (384-well plate) (including blank wells) of the Fatty Acid dye-loading solution.
3. 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.

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

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

EXAMPLE DATA ANALYSIS AND FIGURES

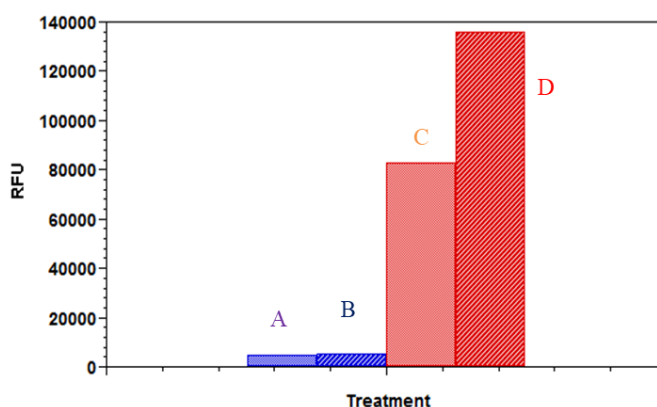


Figure 1. Comparison of fatty acid uptake by 3T3-L1 adipocytes and fibroblast. Cells were plated at 50,000 cells/100 mL/well in a 96 well black wall/clear bottom poly-D lysine plate for 5 hours, and then serum deprived for 1 hour. Cells were treated without (control) or with insulin (150 nM), and incubated at 37 °C, 5% CO₂ 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 FlexStation plate reader using bottom read mode. A – fibroblasts (Control); B – fibroblasts (Insulin); C – adipocytes (Control); D– adipocytes (Insulin).

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.