# ab222944 Fatty Acid Oxidation Complete Assay Kit

For the convenient measurement of Fatty Acid Oxidation (FAO) in live cells.

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

#### 1. Overview

Fatty Acid Oxidation Complete Assay Kit (ab222944) offers a complete solution to measure Fatty Acid Oxidation (FAO) in live cells using conventional fluorescence plate readers. This product combines Fatty Acid Oxidation Assay (ab217602) and Extracellular Oxygen Consumption Assay (ab197243) in one practical kit for single purchase. The protocol booklet is a combination of protocol booklet for ab217602 and the protocol booklet for ab197243.

The assay uses the 18C unsaturated fatty acid Oleate as substrate, and includes two FAO modulators, etomoxir and FCCP. Etomoxir, an inhibitor of the carnitine transporter CPT1, prevents Oleate import and thereby limits the supply of reducing equivalents to the ETC, reducing oxygen consumption in turn. The remaining ETC (electron transport chain) activity is driven by non-long chain FAO. FCCP treatment induces maximal ETC activity by dissipating the mitochondrial membrane potential, while the increased demand for reducing equivalents causes a concomitant increase in the FAO activity. If exogenous long-chain fatty acid is unavailable or import is inhibited, FAO activity will be limited.

Fatty acid oxidation (FAO) is the primary metabolic pathway for degradation of fatty acids. Figure 1 gives an overview of long-chain fatty acid activation, import and oxidation. FAO is an important process in many tissues during periods of glucose deprivation. In organs, such as liver and skeletal muscle, FAO can provide over 75% of cellular ATP while in cardiac tissue it can be responsible for up to 90% of cellular energy requirements. FAO is also now acknowledged as a key factor in cancer metabolism and is also implicated in drug-induced microsteatosis.

# ab217602 Fatty Acid Oxidation Assay

For the convenient measurement of Fatty Acid Oxidation (FAO) in live cells when used in combination with Extracellular Oxygen Consumption Assay (ab197243).

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

# **Table of Contents**

1.	Overview	
2.	Protocol Summary	3
3.	Precautions	4
4.	Storage and Stability	4
5.	Limitations	į
6.	Materials Supplied	į
7.	Materials Required, Not Supplied	(
8.	Technical Hints	7
9.	Reagent Preparation	8
10.	Sample Preparation	9
11.	Assay Controls set up	12
12.	Assay Procedure	14
13.	Typical Assays/Data	17
14.	Notes	20

#### 2. Overview

Fatty Acid Oxidation Assay (ab217602) allows the detection of Fatty Acid Oxidation (FAO) in live cells. This product is designed to be used in combination with our <a href="Extracellular Oxygen Consumption Assay">Extracellular Oxygen Consumption Assay</a> (ab197243).

The assay uses the 18C unsaturated fatty acid Oleate as substrate, and includes two FAO modulators, etomoxir and FCCP. Etomoxir, an inhibitor of the carnitine transporter CPT1, prevents Oleate import and thereby limits the supply of reducing equivalents to the ETC, reducing oxygen consumption in turn. The remaining ETC (electron transport chain) activity is driven by non-long chain FAO. FCCP treatment induces maximal ETC activity by dissipating the mitochondrial membrane potential, while the increased demand for reducing equivalents causes a concomitant increase in the FAO activity. If exogenous long-chain fatty acid is unavailable or import is inhibited, FAO activity will be limited.

Fatty acid oxidation (FAO) is the primary metabolic pathway for degradation of fatty acids. Figure 1 gives an overview of long-chain fatty acid activation, import and oxidation. FAO is an important process in many tissues during periods of glucose deprivation. In organs, such as liver and skeletal muscle, FAO can provide over 75% of cellular ATP while in cardiac tissue it can be responsible for up to 90% of cellular energy requirements. FAO is also now acknowledged as a key factor in cancer metabolism and is also implicated in drug-induced microsteatosis.

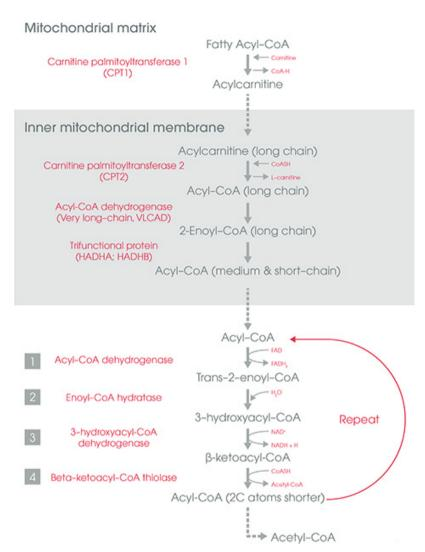


Figure 1. Overview of long-chain fatty acid activation, import and oxidation.

# 3. Protocol Summary

Day 1

Plate cells and return to culture overnight
Prepare Base Measurement Media

Incubate Over Night

Day 2 (OPTIONAL)

Prepare Glucose-Deprivation Media

Replace culture media with Glucose-Deprivation Media

and return cells to culture

Incubate Over Night

Day 3

Prepare FA-Free & FA Measurement Media Prepare FAO kit controls (Etoxomir, FCCP and BSA) Prepare O<sub>2</sub> Consumption reagent (ab197243)



Wash cells twice with FA-Free Measurement Media
Add FA/FA-Free Measurement Media and O<sub>2</sub> consumption reagent
Add controls (Etoxomir, FCCP, BSA)

Overlay plate with High Sensitivity mineral oil (ab197243)



Measure on fluorescence plate reader

Analyze kinetic data output to determine FAO-driven ETC activity

#### 4. Precautions

Please read these instructions carefully prior to beginning the assay.

- All assay kit components have been formulated and quality control tested to function successfully as a kit.
- We understand that, occasionally, experimental protocols might need to be modified to meet unique experimental circumstances.
   However, we cannot guarantee the performance of the product outside the conditions detailed in this protocol booklet.
- Reagents should be treated as possible mutagens and should be handle with care and disposed of properly. Please review the Safety Datasheet (SDS) provided with the product for information on the specific components.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipette by mouth. Do not eat, drink or smoke in the laboratory areas.
- All biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.

# 5. Storage and Stability

Kit has a storage time of 1 year from receipt. Please observe storage conditions of each individual component described in the Materials Supplied section for correct storage upon receipt.

 $\Delta$  Note: Reconstituted reagents are stable for 3 months.

# 6. Limitations

- Kit intended for research use only. Not for use in diagnostic procedures.
- Do not mix or substitute reagents or materials from other kit lots or vendors. Kits are QC tested as a set of components and performance cannot be guaranteed if utilized separately or substituted.

# 7. Materials Supplied

Item	Quantity	Storage temperatur e (before prep)	Storage temperatur e (after prep)
FAO Conjugate (3 mM)	1 mL	4°C	4°C
FAO Control (1.5 mM)	500 µL	4°C	4°C
FAO Tablet (base media)	1 tablet	RT	4°C
L-Carnitine	1 vial	4°C	-20°C
FCCP	1 vial	-20°C	-20°C
Etomoxir	1 vial	-20°C	-20°C

# 8. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- Extracellular Oxygen Consumption Assay (ab197243)
- Microplate reader capable of measuring fluorescence, with suitable filter and plate temperature control – see Instrument and Measurement Settings section on the Extracellular Oxygen Consumption Assay (ab197243) protocol for suitable plate readers
- MilliQ water or other type of double distilled water (ddH<sub>2</sub>O)
- DMSO
- Pipettes and pipette tips, including multi-channel pipette
- Assorted glassware for the preparation of reagents and buffer solutions
- Tubes for the preparation of reagents and buffer solutions
- Sterile 96-well plate (black wall with clear flat bottom), or standard clear plates for cell culture
- Cell culture media
- Base glucose deprivation media: glucose-free DMEM media, 1 mM glucose, 1 mM L-glutamine, 1% FBS, Penicillin/streptomycin solution (100 U/mL /0.1 mg/mL)
- Glucose: to prepare FA-Free Measurement Media
- (Optional) Plate block heater for plate preparation.

To prepare Base measurement media:

- HCl and NaOH: to bring media to pH7.4

#### 9. Technical Hints

- This kit is sold based on number of tests. Number of tests based on described procedure, stock dilutions, concentrations and wash steps. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions.
- Selected components in this kit are supplied in surplus amount to account for additional dilutions, evaporation, or instrumentation settings where higher volumes are required. They should be disposed of in accordance with established safety procedures.
- Avoid foaming or bubbles when mixing or reconstituting components.
- Avoid cross contamination of samples or reagents by changing tips between sample and reagent additions.
- Ensure plates are properly sealed or covered during incubation steps.
- Ensure all reagents and solutions are at the appropriate temperature before starting the assay.
- Make sure all necessary equipment is switched on and set at the appropriate temperature.
- Refer to Instrument and Measurement Settings Table (Table 1) in the user manual for Extracellular O<sub>2</sub> Consumption Assay (ab197243) for recommended settings for your plate reader.

# 10. Reagent Preparation

Briefly centrifuge small vials at low speed prior to opening.

# 10.1 FAO Conjugate (Oleate-BSA conjugate, 3 mM) (1mL):

Ready to use as supplied. Equilibrate to room temperature prior to use. Store at 4°C.

#### 10.2 FAO Control (BSA, 1.5 mM) (500 μL):

Ready to use as supplied. Equilibrate to room temperature prior to use. Store at 4°C.

#### 10.3 FAO Tablet (1 Base media tablet):

To prepare Base Measurement Media, dissolve tablet provided in 100 mL ddH $_2$ O and warm solution to  $37^{\circ}$ C to ensure table is completely dissolved. Adjust pH to 7.4 using HCl and NaOH. Filter sterilize base media. Filtered Base Measurement Media can be stored at  $4^{\circ}$ C for 3 weeks.

#### 10.4 L-Carnitine (L-carnitine hydrochloride, 4 mg):

Prepare a 50 mM L-Carnitine stock solution (100X) by dissolving vial contents in 400  $\mu$ L ddH<sub>2</sub>O. Aliquot L-Carnitine stock solution so that you have enough volume to perform the desired number of assays (recommendation: 100  $\mu$ L to cover 100 tests). Store at -20°C. Once reconstituted, use component within 3 months.

# 10.5 FCCP (0.004 mg):

Prepare a 250  $\mu$ M FCCP stock solution (100X) by dissolving vial contents in 60  $\mu$ L of DMSO. Aliquot FCCP stock solution so that you have enough volume to perform the desired number of assays (recommendation: 20  $\mu$ L to cover 20 tests). Store at - 20°C. Once reconstituted, use component within 3 months.

#### 10.6 Etomoxir (0.074 mg):

Prepare a 400  $\mu$ M Etomoxir stock solution (10X) by dissolving vial contents in 550  $\mu$ L ddH<sub>2</sub>O. Aliquot Etomoxir stock solution so that you have enough volume to perform the desired number of assays (recommendation: 50  $\mu$ L or 100  $\mu$ L to cover 10 or 20 tests respectively). Store at -20°C. Once reconstituted, use component within 3 months.

# 11. Sample Preparation

#### General Sample Information:

- Prepare a cell titration experiment to identify a suitable cell density for a specific cell type and conditions.
- Cells are seeded at a density to achieve full confluence on the day of measurement. Plating density, cell type and basal metabolic rate will determine oxygen consumption rate.
- If performing an overnight glucose deprivation step using nonterminally differentiated cells, seeding densities should be adjusted downwards to facilitate doubling.
- We recommend following the plate map suggested in figure 2 to facilitate assay procedure.

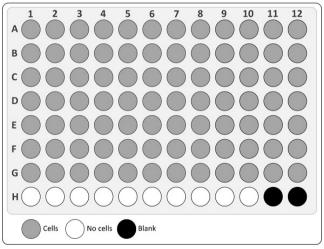


Figure 2. Plate Map.

 If using extended culture periods (> 2 days), we recommend following plate map (figure 3), adding 200 µL of culture medium to all outer wells. This minimizes plate effects related to inconsistent cell growth across the microplate.

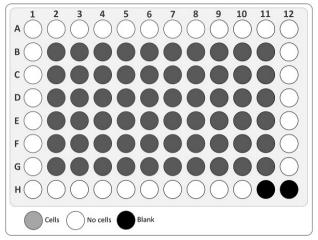


Figure 3. Recommended plate map for extended culture times (> 2 days)

 Inconsistent growth in some cell types can be additionally reduced by allowing the plate to stand at RT for 30 minutes after plating before returning plate to cell culture.

#### 11.1 Adherent cells (2D cell culture):

- 11.1.1 Seed cells in a 96-well plate a density of 6 x  $10^4$  cells/well in 200  $\mu L$  culture medium.
- 11.1.2 Incubate overnight in a CO<sub>2</sub> incubator at 37°C (typical incubation time > 14 hours).
- 11.1.3 OPTIONAL: Glucose deprivation step.

 $\Delta$  Note: Including a glucose deprivation step before performing the assay increases cellular dependence on FAO. For maximum FAO dependence, concentrations of L-Carnitine, glucose and FAO Conjugate should be optimized for each cell type.

- 11.1.3.1 Prepare Base glucose deprivation media as described in Section 7 (media can be stored for 2 weeks at 4°C).
- 11.1.3.2 Prepare Glucose deprivation media by adding 0.5 mM L-Carnitine (1:100 of stock) to Base glucose deprivation media.
- 11.1.3.3 Wash cells twice with Base glucose deprivation media.
- 11.1.3.4 Add 200 µL of Glucose deprivation media.
- 11.1.3.5 Incubate overnight in a  $CO_2$  incubator at 37°C (typical incubation time > 14 hours).

#### 11.2 3D cell culture:

11.2.1 Plate cells for 3D cultures at a higher density than optimized for 2D cultures. When plating 3D cultures, prepare the plate or matrix solution in advance as per manufacturer's instructions.

 $\Delta$  **Note:** test compounds can take longer to penetrate and affect 3D cultures and longer treatment times may be necessary.

11.2.2 OPTIONAL: Glucose deprivation step.

**Δ Note:** Including a glucose deprivation step before performing the assay increases cellular dependence on FAO. For maximum FAO dependence, concentrations of L-Carnitine, glucose and FAO Conjugate should be optimized for each cell type.

- 11.2.2.1 Prepare Base glucose deprivation media as described in Section 7 (media can be stored for 2 weeks at 4°C).
- 11.2.2.2 Prepare Glucose deprivation media by adding 0.5 mM L-Carnitine (1:100 of stock) to Base glucose deprivation media.
- 11.2.2.3 Wash cells twice with Base glucose deprivation media.
- 11.2.2.4 Add 200 µL of Glucose deprivation media.
- 11.2.2.5 Incubate overnight in a  $CO_2$  incubator at 37°C (typical incubation time > 14 hours).

# 12. Assay Controls set up

General guidelines for optimal concentrations and treatment times for assay controls:

#### 12.1 FAO modulators:

- FCCP exhibits a bell-shaped dose-response which can vary between cell types. The concentration which delivers maximum respiratory activity should be titrated for each cell type: run a FCCP serial dilution (15-0.5 µM) in the presence of FAO Conjugate. Higher FCCP concentrations may be required when using FAO Conjugates as compared with glucose-based measurement due to the ability of BSA to bind FCCP (FAO Conjugate is a 2:1 Oleate-BSA conjugate).
- FAO Conjugate is typically used at 150 μM. However, the concentration at which maximum respiratory activity is observed can be cell type dependent. Optimum concentration can be determined by measuring oxygen consumption at varying FAO Conjugate concentrations (typically 50-200 μM) in the presence of FCCP. Users may also wish to add FAO Conjugate to Glucose Deprivation Media (typically 100 μM).
- L-Carnitine is typically used at 0.5 mM. However, the optimum concentration to facilitate LCFA transport is cell type-dependent.
   Optimum concentration can be determined by measuring oxygen consumption at varying L-Carnitine concentrations in the presence of FCCP.
- Etomoxir can exhibit 'off-target' effects if used at > 40 μM. Etomoxir efficacy can be reduced in presence of high serum and BSA concentrations. In these situations, use higher Etomoxir concentrations to ensure CPT-1 inhibition. A minimum of 10 minutes should elapse between Etomoxir treatment and the commencement of measurement to ensure CPT-1 inhibition has impacted oxygen consumption prior to measurement. To maximize inhibition, Etomoxir can be pre-incubated in FA-Free Media prior to the addition of FAO Conjugate or BSA control.

## 12.2 Optional additional controls:

- Antimycin A can be used as optional negative biological control.
   Antimycin blocks the ETC thereby inhibiting ETC-related oxygen consumption. To use in the assay, add 1 μL of Antimycin A 100 μM stock solution (in DMSO) to wells.
- Oligomycin can be used as optional coupling control. Oligomycin blocks the F1/F0 ATPase highlighting the portion of  $O_2$  consumption driving aerobic ATP production. Remaining  $O_2$  consumption is typically due to uncoupled mitochondria. To use in this assay, add 1  $\mu$ L of Oligomycin 100  $\mu$ M stock solution (in DMSO) to wells.

Control	Typical concentration
FCCP	Titration recommended to establish best concentration
FAO-Conjugate	150 μΜ
L-Carnitine	0.5 mM
Etomoxir	< 40 µM
Antimycin A	1 μΜ
Oligomycin	1 μΜ

# 13. Assay Procedure

- This assay is designed to be used as companion kit together in combination with Extracellular Oxygen Consumption Assay (ab197243) The Extracellular Oxygen Consumption Assay User Manual describes instrument set-up, assay optimization, data analysis and troubleshooting. The described instrument set-up and signal optimization steps should be performed prior running an FAO assay.
- We recommend that you assay all controls and samples in triplicate.
- Use a plate block heater for plate preparation and pre-warm plate reader to measurement temperature (typically 37°C).
- Compounds are typically added immediately pre-treatment to determine their effect on FAO and related mitochondrial functional. For some 3D models, a pre-incubation step can be incorporated to ensure compounds access cells within the 3D construct. Longer treatment times can be used as required: in these instances, compound should be present in both culture media (during incubation) and measurement media.
- Long-term measurement with CO<sub>2</sub> control: additional media buffering capacity is required when conducting long term measurements (> 2 h) outside 5% CO<sub>2</sub>. This is achieved by supplementing Measurement Media with 5 mM HEPES.
   Supplementation is not required if plate reader has environmental gas control where 5% CO<sub>2</sub> can be maintained within the measurement chamber.

# 13.1 Prepare additional reagents:

13.1.1 FA-Free Measurement Media: To Base Measurement Media (Step 9.3), add 0.5 mM L-Carnitine (1:100 final dilution) and 2.5 mM glucose.

 $\Delta$  **Note:** Optimal L-Carnitine and glucose concentration may be cell-type specific and maybe require additional optimization.

 $\Delta$  Note: If long term measurements are being performed outside 5% CO<sub>2</sub> a HEPES supplement is recommended.

13.1.2 FA Measurement Media: FA-free Measurement Media + 150 μM FAO Conjugate (1:20 dilution from 3 mM stock – Step 9.1).

13.1.3 Extracellular  $O_2$  Consumption Reagent: prepare reagent as described in Extracellular  $O_2$  Consumption Assay (ab197243) user manual.

#### 13.2 Wash cells:

- 13.2.1 Place the plate on a plate block heater set to assay temperature (typically 37°C) and remove spent culture media with an aspirator (be careful not to dislodge cells from the base of the wells).
- 13.2.2 With a multichannel or repeater pipette, add 100 µL of the prewarmed FA-Free Media to each well.
- 13.2.3 Repeat wash step one more time.

#### 13.3 Add assay media to wells:

- Signal control wells (wells with no cells; row H) = 90 µL of prewarmed FA Measurement Media.
- Blank control wells (H11 and H12) = 90 μL of pre-warmed FA Measurement Media.
- Sample wells = 90 µL of pre-warmed FA Measurement Media.
- FA-Free control wells =85 μL of FA-Free Measurement Media + 5 μL of BSA control.

 $\Delta$  Note: FA-Free Measurement Media is used as a control to measure  $O_2$  consumption without FAO Conjugate. BSA control is added to ensure that the free concentrations of test compounds are consistent between FA and FA-Free conditions. BSA concentration in FA-free control wells should be consistent with the BSA concentration in samples containing FAO Conjugate.

- 13.3.1 Add 10  $\mu$ L of Extracellular O<sub>2</sub> Consumption Reagent (Step 12.1.3) to each sample, FA-Free control and signal control wells. Do not add to blank control wells.
- 13.3.2 Add 10 µL of FA Measurement Media to Blank Control wells.

**\DeltaNote:** If measuring a full 96-well plate, we recommend diluting reconstituted ab197243 stock 1 in 10 in the relevant measurement media and, using a multichannel pipette, to add 100  $\mu$ L to each well. Add 100  $\mu$ L of FA Measurement Media (no ab197243 reagent) to the Blank Control wells.

#### 13.4 Treat cells:

Add relevant treatment or compounds to cells (see Section 13 for more detailed information on how to perform assays). The procedure below describes how to perform drug screening:

13.4.1 Add test compound or vehicle (typically 1-5 µL) to test wells (Step 12.3): we recommend using 6-8 compound dilutions.

 $\Delta$  **Note:** We recommend keeping volume of added compound as low as possible to minimize any potential vehicle effects.

**Δ Note:** Additional BSA control stock is added to wells without Oleate (Step 12.3.1-12.3.2). Cells should be co-treated with FCCP if impact on maximal FAO is being determined. Etomoxir is used as a control.

#### 13.5 Measurement:

- 13.5.1 Seal each well with 100  $\mu$ L of pre-warmed High Sensitivity Mineral Oil (component from ab197243), taking care to avoid bubbles.
- 13.5.2 Read plate immediately in a fluorescence microplate reader as described in the protocol booklet for Extracellular Oxygen Consumption Assay (ab197243).

#### 13.6 Calculations:

13.6.1 Process data as described in Section 15 of the protocol booklet for Extracellular Oxygen Consumption Assay (ab197243).

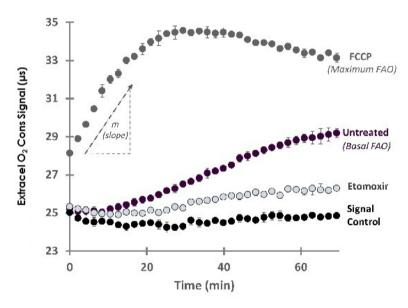
# 14. Typical Assays/Data

#### 14.1 Sample FAO-Driven Oxygen Consumption

Cellular dependence on, or preference for FAO can be determined using Etomoxir and FCCP in the presence or absence of FAO substrate (Oleate). Etomoxir blocks long chain fatty-acid (LCFA) uptake, while FCCP increases cellular energy demand, thereby increasing FAO dependence.

For compound addition, add 10  $\mu$ L of Etomoxir and 1  $\mu$ L FCCP stock to relevant samples on Step 12.3.

 $\Delta$  **Note:** Ensure a minimum of 10 minutes between Etomoxir addition and assay measurement.



**Figure 4.** FAO-driven oxygen respiration in HepG2 cells treated with the CPT-1 inhibitor Etomoxir (white) and uncoupler FCCP (gray).

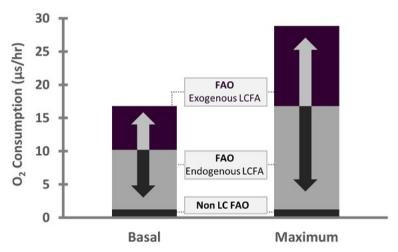
- Untreated cells curve shows a steady increase of the Extracellular
   O<sub>2</sub> Consumption Reagent signal reflecting ETC (electron transfer chain)-driven oxygen consumption.
- **Signal Control** shows probe signal in the absence of cell respiration.

- Etomoxir treatment prevents oleate import, resulting in reduced availability of reducing equivalents and a resultant decrease in ETC activity. The remaining ETC activity (difference between Etoxomir treatment and Signal Control) is driven by metabolic activity other than long chain FAO.
- FCCP treatment induces maximal ETC activity by dissipating the mitochondrial membrane potential. Increased demand for reducing equivalents causes a concomitant increase in FAO as indicated by the rapid increase in Extracellular O<sub>2</sub> Consumption Reagent signal. This strong increase in ETC activity is not observed where exogenous LCFA is unavailable or where import is inhibited.

## 14.2 Evaluating Exogenous and Endogenous FAO

FAO-driven respiratory activity can be investigated further by calculating the rate of signal change for each FAO assay profile, facilitating assessment of exogenous FAO (Oleate supplied), endogenous FAO (Oleate-free) and non-LC FAO (Etomoxir treated). This can be determined using slopes (m) calculated from the linear portion of each profile:

Exogenous FAO =  $m_{Oleate}$  -  $m_{Etomoxir}$ Endogenous FAO =  $m_{Oleate-free}$  -  $m_{Etomoxir}$ Non-LC FAO =  $m_{Etomoxir}$  -  $m_{Signal Control}$ 



**Figure 5:** FAO-driven respiration of HepG2 cells treated with the CPT-1 inhibitor Etomoxir and uncoupler FCCP. The figure summarizes the balance between these parameters under "Basal" and "Maximum" (FCCP treated) conditions. The increased energy demand imposed by FCCP treatment is met by increased FAO.

# 15. Notes

# ab197243 Extracellular O<sub>2</sub> Consumption Assay

For the measurement of extracellular oxygen consumption in isolated mitochondria, cell populations, 3D culture models, tissues and enzymes.

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

# **Table of Contents**

1.	Overview	23
2.	Protocol Summary	24
3.	Precautions	25
4.	Storage and Stability	25
5.	Limitations	26
6.	Materials Supplied	26
7.	Materials Required, Not Supplied	27
8.	Technical Hints	28
9.	Reagent Preparation	29
10.	Plate Reader Set-Up	30
11.	Signal Optimization	36
12.	Sample Preparation	37
13.	Assay Procedure	39
14.	Assay Procedure for 384 well plate	41
15.	Calculations	42
16.	Typical Data	43
17.	Assay Throughput and Performance	44
18.	Additional Assays/Data	46
19.	Notes	51

#### 1. Overview

Extracellular  $O_2$  Consumption Assay (ab197243) measures extracellular oxygen consumption rate (OCR) in a variety of samples.

Mitochondrial dysfunction is implicated in numerous disease states and is also a major mechanism of drug-induced toxicity. Oxygen consumption is one of the most informative and direct measures of mitochondrial function. Traditional methods of measuring oxygen consumption are hampered by the limitations of low throughput and high complexity. The Extracellular  $O_2$  Consumption Assay (ab197243) solves these limitations by providing a direct, real-time measurement of extracellular oxygen consumption rate (OCR) to analyze cellular respiration and mitochondrial function.

The assay is based on the ability of oxygen to quench the excited state of Extracellular  $O_2$  Consumption reagent present in the kit. As the test material respires, oxygen is depleted in the surrounding environment, which is seen as an increase in phosphorescence signal. The addition of a high-sensitivity mineral oil is used to limit back diffusion of ambient oxygen. Measured on standard fluorescence plate readers (96- or 384- well), with standard cell culture microplates, the Extracellular  $O_2$  Consumption Assay (ab197243) is suitable for use with whole cell populations (both adherent and suspension cells), isolated mitochondria, a wide range of 3D culture models, tissues, small organisms, as well as isolated enzymes, bacteria, yeasts and molds.

The flexible plate reader format, allows multiparametric or multiplex combination with other similar products. For example, in combination with Glycolysis Assay (ab197244), the Extracellular  $O_2$  Consumption Assay (ab197243) allows simultaneous real-time measurement of mitochondrial respiration and glycolysis and the analysis of the metabolic phenotype of cells and the shift (flux) between the two pathways under pathological states.

# 2. Protocol Summary

Prepare O<sub>2</sub> Consumption Reagent

Add Reagent (and compounds of interest) to sample

Add high-sensitivity mineral oil

Measure fluorescence in plate reader

#### 3. Precautions

Please read these instructions carefully prior to beginning the assay.

- All kit components have been formulated and quality control tested to function successfully as a kit.
- We understand that, occasionally, experimental protocols might need to be modified to meet unique experimental circumstances. However, we cannot guarantee the performance of the product outside the conditions detailed in this protocol booklet.
- Reagents should be treated as possible mutagens and should be handle with care and disposed of properly. Please review the Safety Datasheet (SDS) provided with the product for information on the specific components.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth. Do not eat, drink or smoke in the laboratory areas.
- All biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.

# 4. Storage and Stability

Store kit at 4°C in the dark immediately upon receipt. Kit has a storage time of 1 year from receipt, providing components have not been reconstituted.

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Materials Supplied section.

Aliquot components in working volumes before storing at the recommended temperature.

 $\Delta$  Note: Reconstituted reagent is stable for 1 month.

# 5. Limitations

- Assay kit intended for research use only. Not for use in diagnostic procedures.
- Do not mix or substitute reagents or materials from other kit lots or vendors. Kits are QC tested as a set of components and performance cannot be guaranteed if utilized separately or substituted.

# 6. Materials Supplied

Item	Quantity	Storage condition (before prep)	Storage condition (after prep)
Extracellular O <sub>2</sub> Consumption Reagent	1 vial	4°C	-20°C
High Sensitivity Mineral Oil	1 dropper bottle	4°C	RT

# 7. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- Microplate reader capable of measuring fluorescence, with suitable filter and plate temperature control – see Instrument and Measurement Settings section for suitable plate readers
- Double distilled water (ddH<sub>2</sub>O)
- Pipettes and pipette tips, including multi-channel pipette
- Assorted glassware for the preparation of reagents and buffer solutions
- Tubes for the preparation of reagents and buffer solutions
- Sterile 96 well plate (black wall with clear flat bottom), or standard clear plates for cell culture
   For cells:
- Cell culture medium
   For isolated mitochondria:
- Measurement buffer (250 mM sucrose, 15 mM KCl, 1 mM EGTA, 5 mM MgCl $_2$ , 30 mM, K $_2$ HPO $_4$ , pH 7.4)
- Mitochondrial substrate (succinate, glutamate or malate)
- ADP

#### 8. Technical Hints

- This kit is sold based on number of tests. A "test" simply refers to a single assay well. The number of wells that contain sample, control or standard will vary by product. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions.
- Selected components in this kit are supplied in surplus amount to account for additional dilutions, evaporation, or instrumentation settings where higher volumes are required. They should be disposed of in accordance with established safety procedures.
- Avoid foaming or bubbles when mixing or reconstituting components.
- Avoid cross contamination of samples or reagents by changing tips between sample, standard and reagent additions.
- Ensure plates are properly sealed or covered during incubation steps.
- Ensure all reagents and solutions are at the appropriate temperature before starting the assay.
- Make sure all necessary equipment is switched on and set at the appropriate temperature.
- Refer to Instrument and Measurement Settings table (Table 1) for recommended settings for your plate reader.
- While compatible with all plate types, black border clear bottom plats give optimal signal-to-noise ratios.
- For first time users, we recommend performing a Signal Optimization Step (see Section 11).

# 9. Reagent Preparation

Briefly centrifuge small vials at low speed prior to opening.

#### 9.1 Extracellular O<sub>2</sub> Consumption Reagent:

Prepare a stock solution of the Extracellular  $O_2$  Consumption Reagent by adding 1 mL of ddH<sub>2</sub>O, PBS, culture media or buffer to the vial. Mix by gently aspirating 3-4 times.

Recommended working dilution = 1/15 (10 µL per 150 µL of sample for 1x 96-wp).

Avoid freeze/thaw. Reconstituted reagent is stable for one month.

#### 9.2 High Sensitivity Mineral Oil:

Ready to use as supplied. Pre-warm to 37°C prior to use.

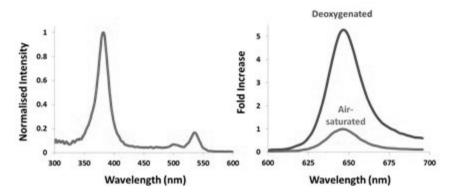
Although mineral oil is provided in a dropper bottle for convenience, we recommend using a repeater pipette for routine use. To apply oil using a repeater pipette, trim 3 – 4 mm off the tip at a 45° angle. Remove internal nozzle cap from the dropper bottle and slowly pick up the pre-warmed mineral oil.

Store mineral oil at room temperature in the dark.

# 10. Plate Reader Set-Up

#### 10.1 Measurement Parameters

The Extracellular Consumption Reagent is a chemically stable and inert, biopolymer-based, cell impermeable oxygen-sensing fluorophore.



**Figure 1.** Excitation and emission spectra of the Extracellular  $O_2$  Consumption Reagent. Left panel shows normalized excitation (Ex = 360 – 400 nm; Peak 380 nm). Right panel shows emission (Em = 630 – 680 nm; Peak 650 nm) in deoxygenated and oxygenated (air-saturated) conditions.

	Peak Maxima (nm)	Peak (nm)
Excitation*	380	360 – 400
Emission	650	630 – 680

<sup>\*</sup>Excitation at  $532 \pm 7.5$  nm is also possible

- This assay is measured with prompt or time-resolved fluorescence (TR-F) readers, monochromator or filter-based. Optimal wavelengths are  $\lambda = 380$  nm for excitation ( $\lambda = 532$  nm can also be used) and  $\lambda = 650$  nm for emission.
- Probe signals should be at least 3 times above blank signal.
- The O<sub>2</sub> probe response is temperature dependent, so good temperature control of the plate during the measurement is important.

#### 10.2 Fluorescence measurements

Outlined below are three fluorescence modalities that can be used with this assay, depending on the plate reader type and instrument setup.

10.2.1 Basic: Intensity Measurement

Measurement of Signal Intensity (sometimes referred to as Prompt) provides flexibility to use wide range of commonly available fluorescence-, monochromator or filter-based plate readers. Optimal wavelengths are  $\lambda$  = 380 nm for excitation and  $\lambda$  = 650 nm for emission, with detection gain parameters (PMT) typically set at medium or high.

**\Delta Note:** Extracellular consumption reagent should return Signal to Blank ratio (S:B)  $\geq$  3.

#### 10.2.2Standard: TR-F Measurement

Using time-resolved fluorescence (TR-F) will increase performance levels. TR-F measurement reduces non-specific background and increases sensitivity.

Optimal delay time is 30 µs and gate (integration) time is 100 µs.

**\Delta Note:** Extracellular consumption reagent should return Signal to Blank ratio (S:B)  $\geq$  3. S:B  $\sim$  10 are typical.

## 10.2.3 Advanced: Dual-Read TR-F (Lifetime calculation)

Optimal performance can be achieved using dual-read TR-F in combination with subsequent ratiometric Lifetime calculation, to maximize dynamic range.

**\Delta Note:** Extracellular consumption reagent should return Signal to Blank ratio (S:B)  $\geq$  3. S:B up to 60 are possible.

Dual-read TR-F and subsequent Lifetime calculation allows measurement of the rate of fluorescence decay of the Extracellular consumption reagent, and can provide measurements of oxygen consumption that are more stable and with a wider dynamic range than measuring signal intensity.

Optimal dual-delay and gate (integration) times:

- Integration window 1: 30 μs delay (D1), 30 μs measurement time (W1)
- Integration window 2: 70 μs delay (D2), 30 μs measurement time (W2)

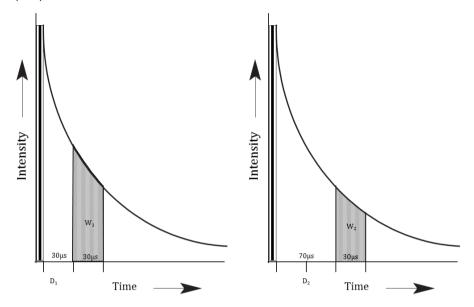


Figure 2. Illustrating dual read TR-F measurement.

Use the dual intensity readings to calculate the corresponding Lifetime (µs) using the following transformation:

Lifetime ( $\mu$ s) [T] = (D2-D1)/ln(W1/W2)

Where W1 and W2 represent the two (dual) measurement windows and D1 and D2 represent the delay time prior to measurement of W1 and W2 respectively. This provides Lifetime values in µs at each measured time point for each individual sample (Figure 2).

**\Delta Note:** S:B for Integration window 2 is recommended to be  $\geq$  10 to allow accurate Lifetime calculation. Range of Lifetime values should be 22 – 68 µs, and should only be calculated from samples containing reagent. Lifetime values should not be calculated from blank wells.

See Instrument and Measurement Setting table below for instrument-specific setting and filters. Readers equipped with a TR-F mode, may achieve improved performance using delay and gate time of 30  $\mu$ s and 100  $\mu$ s.

**Table 1.** Recommended Instrument and Measurement Settings

Instrument	Optical Configuration	Intg1 (D <sub>1</sub> /W <sub>1</sub> )	Optimum Mode	Ex (nm)
		Intg2 (D <sub>2</sub> /W <sub>2</sub> )		Em (nm)
BioTek:	Filter-based	30 / 30µs	Dual-read TR-F	Ex 380 ± 20nm Em 645 ± 15nm
Cytation 3 / 5	Top or bottom read	70 / 30µs	(Lifetime)	
BioTek:	Filter-based	30 / 30µs	Dual-read TR-F	Ex 380 ± 20nm
Synergy H1, Neo, 2	Top or bottom read	70 / 30µs	(Lifetime)	Em 645 ± 15nm
BMG Labtech:	Filter-based	30 / 30µs	Dual-read TR-F	Ex 340 ± 50nm (TR-EX)
CLARIOstar	Bottom read	70 / 30μs	(Lifetime)	Em 665 ± 50nm or Em 645± 10nm With LP-TR Dichroic
BMG Labtech:	Filter-based	30 / 30µs	Dual-read TR-F	Ex 340 ± 50nm (TR-EXL)
FLUOstar Omega /	Top or bottom read	70 / 30µs	(Lifetime)	Em 655 ± 25nm (BP-655)
POLARstar Omega				
Perkin Elmer:	Filter-based	30 / 30µs	Dual-read TR-F	Ex 340 ± 40nm (D340)
VICTOR series/ X4, X5	Top read	70 / 30µs	(Lifetime)	Em 642 ± 10nm (D642)

Instrument	Optical Configuration	Intg1 (D <sub>1</sub> /W <sub>1</sub> ) Intg2 (D <sub>2</sub> /W <sub>2</sub> )	Optimum Mode	Ex (nm) Em (nm)
		9_ (-2,2)		
Tecan:	Monochromator /	30 / 30µs	Dual-read TR-F	Ex 380 ± 20nm
Infinite M1000Pro /	Filter-based	70 / 30µs	(Lifetime)	Em 650 ±20nm or Em 670±40nm
F200Pro	Top or bottom read			
BioTek:	Monochromator /	30 / 100µs	TR-F	Ex 380±20nm
Synergy HTx / Mx	Filter-based	n/a		Em 650±15nm
	Top or bottom read			
BMG Labtech:	Filter-based	40 / 100µs	TR-F	Ex 337 nm (HTRF Module)
PHERAstar FS	Top or bottom read	n/a		Em 665 nm (HTRF Module)
BMG Labtech:	Filter-based	30 / 100µs	TR-F	Ex 340 ± 50nm (TR-EXL)
FLUOstar Optima /	Top or bottom read	n/a		Em 655 ± 50nm (BP-655)
POLARstar Optima				

Instrument	Optical Configuration	Intg1 (D <sub>1</sub> /W <sub>1</sub> ) Intg2 (D <sub>2</sub> /W <sub>2</sub> )	Optimum Mode	Ex (nm) Em (nm)
Perkin Elmer:	Filter-based	40 / 100µs	TR-F	Ex 340 ±60nm (X340)
EnVision	Top read	n/a		Em 650 ± 8nm (M650)
Perkin Elmer:	Monochromator based	40 / 100µs	TR-F	Ex 380 ±20nm
EnSpire	Top read	n/a		Em 650±20nm
Tecan:	Monochromator /	30 / 100µs	TR-F	Ex 380±20nm
Infinite M200Pro / Saffire / Genios Pro	Filter-based	n/a		Em 650±20nm
Genios Pro	Top or bottom read			
Mol. Devices:	Monochromator based	n/a	Intensity (Prompt)	Ex 380nm
SpectraMax / Flexstation / Gemini	Top or bottom read	n/a		Em 650nm

# 11. Signal Optimization

- This step is recommended for first time users.
- Use a plate block heater for plate preparation and pre-warm plate reader to measurement temperature.
- 11.1 Prepare 8 replicate wells of a 96-well plate, by adding 150 µL pre-warmed culture medium to each well (A1-A4, B1-B4).
- 11.2 Add 10  $\mu$ L reconstituted Extracellular O<sub>2</sub> Consumption Reagent to 4 of the replicate wells (A1-A4) and 10  $\mu$ L water, PBS or media to the remaining replicates wells (B1-B4).
- 11.3 Promptly add two drops (or  $100 \, \mu L$ ) pre-warmed High sensitivity mineral oil to all eight replicate wells, taking care to avoid air bubbles.
- 11.4 Read plate immediately in a fluorescence plate reader over 30 minutes (read every 2-3 minutes.
- 11.5 Examine Signal Control well (A1-A4) and Blank Control well (B1-B4) readings (linear phase) and calculate Signal to Blank (S:B) ratio.

 $\Delta$  Note: For dual read TR-F, calculate S:B for each measurement window.

	1	2	3	4
A	Media + O <sub>2</sub> Reagent + Oil			
В	Media + Oil	Media + Oil	Media + Oil	Media + Oil

# 12. Sample Preparation

### General Sample Information:

- Prepare a cell titration experiment to identify a suitable cell density for a specific cell type and conditions.
- Prepare test compounds for sample treatment as desired.
   Example of typical compounds that can be used as assay control are shown in the table below.

Typical control	Stock concentration
Antimycin A (Complex III inhibitor)	150 μM in DMSO
FCCP (ETC uncoupler)	Titration recommended to establish best concentration
Glucose Oxidase (positive signal control)	1 mg/mL in ddH <sub>2</sub> O

#### 12.1 Adherent cells:

- 12.1.1 Seed cells in a 96-well plate at a density of  $4-8 \times 10^5$  cells/well in 200 µL culture medium.
- 12.1.2Incubate overnight in a CO<sub>2</sub> incubator at 37°C.

 $\Delta$  **Note:** Prepare a cell titration experiment to identify a suitable cell density for a specific cell type and conditions.

## 12.2 Suspension cells:

Seed cells in a 96-well plate at a density of 4 x  $10^6$  cells/well in  $150~\mu L$  culture medium.

 $\Delta$  **Note:** Prepare a cell titration experiment to identify a suitable cell density for a specific cell type and conditions.

#### 12.3 Isolated mitochondria:

 $\Delta$  **Note:** Mitochondria should be freshly prepared as per user's protocol and should not be left on ice longer than recommended in the literature.

Prepare a cell titration experiment to identify a suitable cell density for a specific cell type and conditions.

Initial isolated mitochondria assay optimization: prepare a six-point dilution series of mitochondrial preparation in respiration buffer in 1.5 mL total volume for each concentration.

- 12.3.1 Prepare measurement buffer as follows: 250 mM sucrose, 15 mM KCl, 1 mM EGTA, 5 mM MgCl<sub>2</sub>, 30 mM, K<sub>2</sub>HPO<sub>4</sub>; adjust to pH 7.4.
- 12.3.2Dilute isolated mitochondria to the desired concentration (typical range = 0.125 1.5 mg/mL final concentration) in measurement buffer, depending on the substrate(s) used and which respiration state is being measured.

# 13. Assay Procedure

- We recommend that you assay all controls and samples in duplicate.
- Prepare all controls and samples as directed in the previous sections.
- Use a plate block heater for plate preparation and pre-warm plate reader to measurement temperature (typically 37°C; 30°C for mitochondria).
- Sufficient cell numbers are required to produce measurable signal changes. Oxygen consumption rate is cell-type dependent – highly glycolytic cells may need to be trypsinized and concentrated prior to measurement.

#### PROTOCOL FOR CELLS:

## 13.1 Plate loading:

- 13.1.1 Adherent cells: remove culture media from all assay wells and replace with 150 µL of fresh culture media.

  Suspension cells: ready to use as prepared in Step 12.2.
- 13.1.2Blank controls (we suggest using wells H11 and H12): add 150  $\,$  µL fresh culture media.

## 13.2 Assay set up:

- 13.2.1 Add 10  $\mu$ L reconstituted Extracellular  $O_2$  Consumption Reagent to each sample well.
- 13.2.2Add 10 µL of fresh culture media to blank control wells.
- 13.2.3Add 1 10  $\mu$ L test compound (vehicle control and/or stock) to the wells.

 $\Delta$  **Note:** we recommend keeping the volume of added compound low to minimize any potential effects of solvent vehicle.

13.2.4Promptly seal each well by adding 100  $\mu$ L (or 2 drops) of prewarmed High Sensitivity mineral oil, taking care to avoid air bubbles.

 $\Delta$  **Note:** plate preparation time should be kept to a minimum.

#### 13.3 Measurement:

- 13.3.1 Insert the prepared plate into a fluorescence plate reader pre-set to the measurement temperature (typically 37°C).
- 13.3.2 Measure Extracellular  $O_2$  Consumption signal at 1.5 min intervals for 90 120 minutes (longer for more glycolytic cells) at Ex/Em = 380/650 nm.

#### PROTOCOL FOR ISOLATED MITOCHONDRIA:

- 13.4 Dilute reconstituted Extracellular  $O_2$  Consumption Reagent (Step 9.1)1:10 in measurement buffer.
- 13.5 Add 100 µL reconstituted probe to each sample well.
- 13.6 Add 1 µL test compound in appropriate solvent to the wells.
- 13.7 Add 50 µL of diluted isolated mitochondria (Step 12.3) to each test well. For blank control wells (wells H11 and H12), add 200 µL fresh culture media.
- 13.8 Dissolve substrate in measurement buffer and add 50  $\mu$ L of solution to test wells (see table below for suggested concentrations). Do not add substrate to blank control wells.

Substrate	Mitochondria Concentration (mg/mL)	Typical final substrate concentration (mM)			
Basal state [State 2]					
Glutamate/Malate	1.5	12.5 / 12.5			
Succinate	1.0	25			
ADP-stimulated respiration rate [State 3]					
Glutamate/Malate/ADP	1.0	12.5 / 12.5 / 1.65			
Succinate/ADP	0.5	25 / 1.65			

- 13.9 Promptly seal each well by adding 100 µL (or 2 drops) of High Sensitivity mineral oil, pre-warmed at 30°C, taking care to avoid bubbles.
- **13.10** Insert the prepared plate into a fluorescence plate reader pre-set to 30°C.
- **13.11** Measure Extracellular  $O_2$  Consumption signal at 1.5 min intervals for 10 30 minutes at Ex/Em = 380/650 nm.

# 14. Assay Procedure for 384 well plate

This kit provides enough reagent to perform 200 tests in 384-wp format (half plate).

Follow the same recommendations outlined in Section 13.

Adherent cells: seed cells in a 384-wp at a density of 2 – 4 x 10<sup>5</sup> cells/well in 100 µL culture medium overnight.

### 14.1 Cell preparation:

- Adherent cells: seed cells in a 384-wp at a density of 2 4 x 10<sup>5</sup> cells/well in 100 µL culture medium overnight.
- Suspension cells: prepare a cell concentration stock of 4 x 10<sup>6</sup> cells/mL. Add 75 µL cells per well.

 $\Delta$  **Note:** Prepare a cell titration experiment to identify a suitable cell density for a specific cell type and conditions.

### 14.2 Plate loading:

- 14.2.1 Adherent cells: remove culture media from all assay wells and replace with 75 µL of fresh culture media.

  Suspension cells: ready to use as prepared in Step 14.1.
- 14.2.2Blank controls (we suggest using wells H11 and H12): add 75 µL fresh culture media (for cell-based assay).

## 14.3 Assay set up:

- 14.3.1 Add 5  $\mu$ L reconstituted Extracellular  $O_2$  Consumption Reagent to each sample well.
- $14.3.2 \text{Add} 5 \, \mu\text{L}$  of fresh culture media to blank control wells.
- 14.3.3Add 1 5  $\mu$ L test compound (vehicle control and/or stock) to the wells.

 $\Delta$  **Note:** we recommend keeping the volume of added compound low to minimize any potential effects of solvent vehicle.

14.3.4Promptly seal each well by adding 50 µL (or 1 drops) of prewarmed High Sensitivity mineral oil, taking care to avoid air bubbles.

 $\Delta$  **Note:** High Sensitivity mineral oil is very viscous and it might be difficult to plate into smaller wells.

 $\Delta$  **Note:** plate preparation time should be kept to a minimum.

#### 14.4 Measurement:

14.4.1 Follow instructions described in page 18.

### 15. Calculations

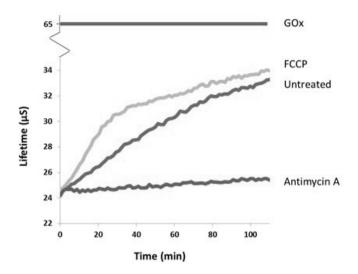
- 15.1 Plot the Blank control well-corrected Extracellular O<sub>2</sub> consumption assay Intensity or Lifetime values versus Time (min).
- 15.2 Select the linear proportion of the signal profile (avoiding any initial lag of subsequent plateau) and apply linear regression to determine the slope (OCR) and correlation coefficient for each well.

 $\Delta$  **Note:** this approach is preferable to calculating a slope from averaged profiles.

15.3 Tabulate the slope values for each test sample, calculating appropriate average and standard deviation values across replicate wells. If optional Signal Control wells are included, the slope obtained for the Signal Control (sample without cells) should be subtracted from all test values.

# 16. Typical Data

Data provided for demonstration purposes only.

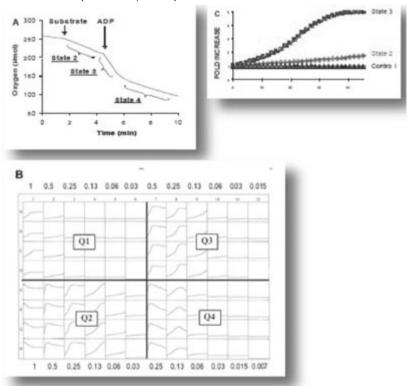


**Figure 3.** Typical Lifetime profile of Extracellular  $O_2$  Consumption Assay for adherent cells, treated with different electron transport chain (ETC) modulator compounds, including antimycin A (recommended negative control). Effect of glucose oxidase (GOx) as positive signal control is illustrated schematically.

# 17. Assay Throughput and Performance

Data output from the Extracellular  $O_2$  Consumption Assay (96 samples) was compared to a polarographic analysis (one sample) of mitochondrial oxygen consumption in Figure 4.

Figure 4A shows typical polarographic analysis illustrating initiation of State 2 (basal) and State 3 (stimulated state) respiration through addition of substrate and ADP respectively. Figure 4B shows oxygen consumption measured using the Extracellular  $O_2$  Consumption Assay, where glutamate/malate- (left panel) and succinate- (right panel) driven oxygen consumption is measured at decreasing mitochondrial protein concentrations in both State 2 (top panel) and State 3 (bottom panel).



**Figure 4.** Analysis of isolated mitochondria using conventional polarograhy (A) and Extracellular  $O_2$  Consumption Assay (n = 4, %CV < 3%) (B). Figure 4C illustrates activation of succinate-driven mitochondrial oxygen consumption measured in figure 4B.

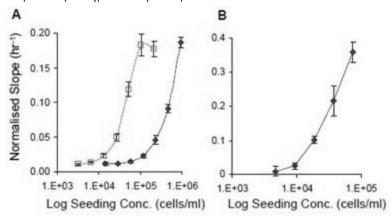
The compatibility of this assay with the microplate format allows analysis under 96 or 384 discrete conditions. The effectiveness of this level of throughput in analyzing isolated mitochondria is highlighted in Figure 4B, which examines increasing mitochondrial protein concentrations on glutamate/malate- and succinate-driven respiration in both basal (State 2) and ADP activated (State 3) states, all tested in quadruplicates.

The performance of the assay is highlighted in Figure 4C, with a coefficient of variance (%CV) below 3%.

# 18. Additional Assays/Data

## 18.1 Monitoring cell respiration

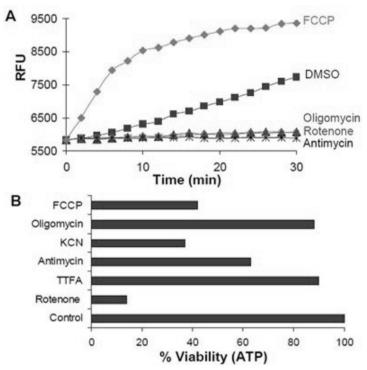
The ability of this product to assess cellular respiration is illustrated in Figure 4. Dilutions curves for HepG2 cells (panel A) and primary rat hepatocytes (panel B) are presented.



**Figure 5.** Cell dilutions measured on 96-well plates using Extracellular  $O_2$  Assay. HepG2 cells (A) were assayed after an overnight (open squares) or 2-day culture period (dark squares); primary rat hepatocytes (B) were assayed after an overnight culture period. Rates of probe signal change (slope of fluorescence signal) were normalized against initial intensity.

## 18.2 Extracellular O<sub>2</sub> Consumption Assay and ATP analysis of cells

To contrast the sensitivities of oxygen consumption and cellular ATP concentrations as indicators of mitochondrial dysfunction, HepG2 cells were treated with a panel of classical ETC modulators. Both cellular oxygen consumption and cellular ATP concentrations were measured.

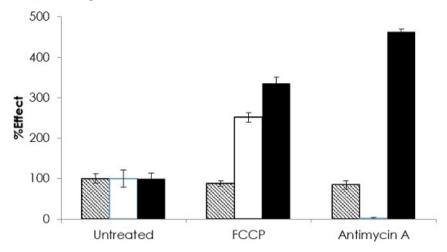


**Figure 6.** Parallel analysis of classical ETC inhibitors on HepG2 cells. 80,000 cells/well were plated and allowed to adhere overnight before performing the assay. Oxygen consumption was measured immediately after treatment, whereas ATP concentration was measured 24 hr post-treatment. Extracellular O<sub>2</sub> consumption data (panel A) illustrates that drug induced mitochondrial dysfunction is evident immediately post-treatment with both inhibition (induced by oligomycin, rotenone or antimycin) and uncoupling (induced by FCCP) being detected. Despite this dysfunction and an additional 24 hr exposure, analysis of cellular ATP concentration (panel B) showed high levels of viability in most scenarios.

This pattern is also replicated in other cell lines and when using other assays.

## 18.3 Cellular energy flux analysis

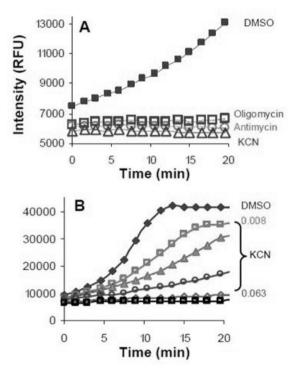
Multiparametric (or multiplex) combination of Extracellular  $O_2$  Consumption Assay together with Glycolysis Assay (ab197244) allows the simultaneous real-time measurement of mitochondrial respiration and glycolysis, leading to the analysis of the metabolic phenotype of cells and the shift (flux) between the two pathways under pathological states.



**Figure 7.** Cellular Energy Flux for HepG2 cells (seeded at 65,000 per well), treated with a combination of drug compounds modulating the ETC (Antimycin A [1  $\mu$ M] and FCCP [2.5  $\mu$ M]), shown as a percentage relative to untreated control cells. Comparative measurements were taken with Extracellular Oxygen Consumption Assay (ab197243) (white column) and Glycolysis Assay [Extracellular acidification] (ab197244) (black column) show the shift between mitochondrial respiration and glycolysis and the cellular control of energy (ATP; measured 1h post-treatment using Luminescent ATP Detection Assay kit (ab113849) (striped column)).

#### 18.4 Assessment of classical mitochondrial effects

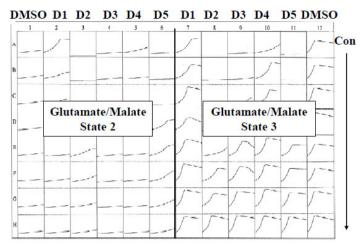
Validation of Extracellular  $O_2$  Consumption Assay for assessment of mitochondria is illustrated in the figure below. The data illustrate the inhibition of mitochondrial function using a panel of classical mitochondrial inhibitors and highlight the dose dependence of this inhibition for KCN (potassium cyanide).



**Figure 8.** Effect of a panel of classical ETC inhibitors on mitochondrial function (A) and dose response inhibition of mitochondrial function by KCN (B).

## 18.5 Compound Screening

Extracellular  $O_2$  Consumption Assay allows screening of compounds at multiple concentrations and in multiple conditions in a single microtiter plate as illustrated in Figure 9. Such data can be processed further to generate dose response data.



**Figure 9.** Effect of test compounds (D01 – D05), on both State 2 and State 3 isolated mitochondrial function using Extracellular  $O_2$  Consumption Assay. Some compounds uncouple in a dose dependent manner while other inhibit.

A 96-well plate format allows screening of 200 compounds a day at a single dose, or acquisition of dose response characteristics for 25 compounds per day. This capability represents a fundamental shift in the capacity for mitochondrial toxicity testing in drug discovery programs, without compromising data quality or information content.

# 19. Notes

# **Technical Support**

Copyright © 2017 Abcam, All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

#### **Austria**

wissenschaftlicherdienst@abcam.com | 019-288-259

#### France

supportscientifique@abcam.com | 01.46.94.62.96

### Germany

wissenschaftlicherdienst@abcam.com | 030-896-779-154

### Spain

soportecientifico@abcam.com | 91-114-65-60

#### **Switzerland**

technical@abcam.com

Deutsch: 043-501-64-24 | Français: 061-500-05-30

UK, EU and ROW

technical@abcam.com | +44(0)1223-696000

#### Canada

ca.technical@abcam.com | 877-749-8807

## US and Latin America

us.technical@abcam.com | 888-772-2226

### Asia Pacific

hk.technical@abcam.com | (852) 2603-6823

#### China

cn.technical@abcam.com | 400 921 0189 / +86 21 2070 0500

## Japan

technical@abcam.co.jp | +81-(0)3-6231-0940

## Singapore

sg.technical@abcam.com | 800 188-5244

#### **Australia**

au.technical@abcam.com | +61-(0)3-8652-1450

### New Zealand

nz.technical@abcam.com | +64-(0)9-909-7829