

Version 2 Last updated 6 January 2020

ab222946 Glycolysis Stress Test Complete Assay Kit

For the characterization of glycolytic stress in live cells.

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

1. Overview

Glycolysis Stress Test Complete Assay Kit (ab222946) offers a complete solution to measure glycolytic characterization at user-defined glucose concentrations in live cells using conventional fluorescence plate readers. This product combines Glycolysis Assay [Extracellular acidification] (ab197244) and Glycolysis Stress Test Companion Assay (ab222945) in one practical kit for single purchase.

The assay uses 2-Deoxyglucose (2-DG) and oligomycin to facilitate an in-depth analysis of cellular glycolytic flux: 2-DG blocks glycolysis through competitive hexokinase inhibition allowing quantification of non-glycolytic extracellular acidification (ECA); oligomycin inhibits ATP synthase, therefore preventing aerobic ATP generation and forcing cells to increase glycolytic flux to meet ATP demand. The levels of such compensatory glycolysis achievable under test conditions can reveal underlying glycolytic perturbation not evident under basal conditions and facilitates assessment of ATP-demand limited glycolysis capability. Glucose oxidase is included in the kit to provide a convenient cell-free ECA assay signal control.

Glycolysis is a critical ATP generating pathway in eukaryotic cells and plays a central role in numerous pathologies including cancer, diabetes and neurodegenerative disorders. The ability to conveniently characterize the glycolytic profile of in vitro cell models used in these disease areas and to determine how specific interventions such as nutrition deprivation, hypoxia or drug treatment impact the profile is therefore of considerable utility.

Version 1 Last updated 6 January 2020

ab222945

Glycolysis Stress Test Companion Assay

For the characterization of glycolytic stress in live cells when used in combination with Glycolysis Assay [Extracellular acidification] (ab197244).

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

Table of Contents

1. Overview	4
2. Protocol Summary	6
3. Precautions	7
4. Storage and Stability	7
5. Limitations	8
6. Materials Supplied	8
7. Materials Required, Not Supplied	9
8. Technical Hints	10
9. Reagent Preparation	11
10. Sample Preparation	12
11. Assay Procedure	14
12. Typical Assays/Data	17
13. FAQs	21
14. Notes	22

1. Overview

Glycolysis Stress Test Companion Assay (ab222945) allows glycolytic characterization at user-defined glucose concentrations in live cells when used in combination with our Glycolysis Assay [Extracellular Acidification] (ab197244).

The assay uses 2-Deoxyglucose (2-DG) and oligomycin to facilitate an in-depth analysis of cellular glycolytic flux: 2-DG blocks glycolysis through competitive hexokinase inhibition allowing quantification of non-glycolytic extracellular acidification (ECA); oligomycin inhibits ATP synthase, therefore preventing aerobic ATP generation and forcing cells to increase glycolytic flux to meet ATP demand. The levels of such compensatory glycolysis achievable under test conditions can reveal underlying glycolytic perturbation not evident under basal conditions and facilitates assessment of **ATP-demand limited glycolysis capability**. Glucose oxidase is included in the kit to provide a convenient cell-free ECA assay signal control.

Glycolysis is a critical ATP generating pathway in eukaryotic cells and plays a central role in numerous pathologies including cancer, diabetes and neurodegenerative disorders. The ability to conveniently characterize the glycolytic profile of *in vitro* cell models used in these disease areas and to determine how specific interventions such as nutrition deprivation, hypoxia or drug treatment impact the profile is therefore of considerable utility.

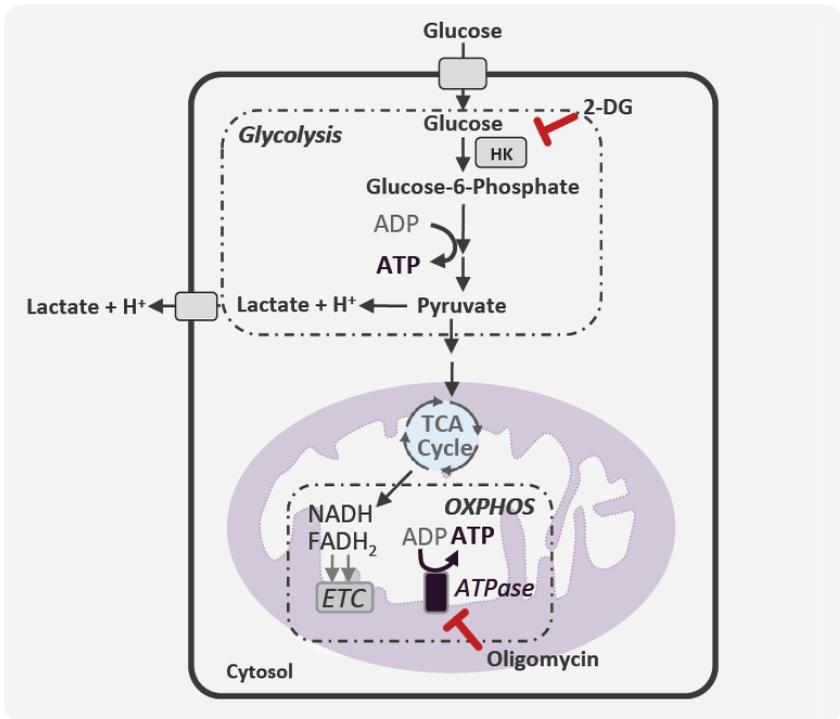


Figure 1. Schematic overview of glycolysis, oxidative phosphorylation (OXPHOS) and the specific targets of the inhibitors provided in the kit.

Glucose is taken up through specific transporters and subsequently converted to pyruvate through a series of reactions, yielding two equivalents of ATP per molecule of glucose use. Pyruvate is then reduced to lactate, eventually leading to the net production of a hydrogen ion (glycolysis). The hydrogen ions generated are exported from the cytoplasm to maintain intracellular pH, in turn, causing extracellular acidification. Pyruvate can also enter the TCA cycle, generating reducing equivalents that are used in the electron transport chain (ETC) to generate a proton gradient across the inner mitochondrial membrane, which in turn drives the generation of ATP catalyzed by F₁/F₀ ATP-synthase (OXPHOS).

2-Deoxyglucose (2-DG) is a competitive inhibitor of hexokinases (HK), and, by extension of, glucose-drive extracellular acidification. Oligomycin inhibits the proton flux through the F₁/F₀ ATP synthase, thereby inhibiting ATP generation via OXPHOS.

2. Protocol Summary

Day 1

Plate cells and return to culture overnight



Incubate Over Night

Day 2

Prepare Glucose-Free Respiration Buffer

Prepare Glycolysis Stress Test controls

Prepare Glycolysis Assay Reagent (ab197244)



Place cells in CO₂-free incubation for 2 hours

Wash cells twice with Glucose-free Respiration Buffer

Add Glucose-free and/or Glucose-containing Respiration Buffer
and Glycolysis Assay Reagent (ab197244)

Add control compounds at pre-optimized concentrations



Measure on fluorescence plate reader using recommended settings

Analyze kinetic data output to determine sample glycolytic
characterization

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 handled 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 -20°C (Respiration Buffer tablet and glucose should be stored at RT) in the dark immediately upon receipt. Kit has a storage time of 1 year from receipt.

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.

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 temperature (before prep)	Storage temperature (after prep)
2-Deoxyglucose (2-DG)	600 µL	-20°C	-20°C
Glucose (720 mg)	1 vial	RT	4°C
Glucose Oxidase (112.5 µg)	1 vial	-20°C	-20°C
Glucose-free Respiration Buffer	1 tablet	RT	4°C
Oligomycin (10.4 µg)	1 vial	-20°C	-20°C
Glycolysis Assay Reagent (lyophilized)	1 vial	4°C	4°C
Respiration Buffer	1 tablet	RT	4°C

7. Materials Required, Not Supplied

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

- Glycolysis Assay [Extracellular acidification] (ab197244)
- Microplate reader capable of measuring fluorescence, with suitable filter and plate temperature control – see Instrument and Measurement Settings section on the Glycolysis Assay [Extracellular acidification] Assay (ab197244) protocol for suitable plate readers
- CO₂-free incubator
- Double distilled water (ddH₂O)
- 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
- Sterilizing filter 0.22 µm
- HCl and KOH – for adjusting pH
- Cell culture medium

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 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.

9. Reagent Preparation

Briefly centrifuge small vials at low speed prior to opening.

9.1 2-Deoxyglucose (2-DG, 500 mM):

Ready to use as supplied. Equilibrate to room temperature before use. Store at -20°C.

9.2 Glucose (720 mg D-(+)-Glucose):

Reconstitute Glucose powder in 4 mL of ddH₂O to generate a 1 mM Glucose stock solution. Filter sterilize using a 0.22 µm filter. Store at 4°C. Use Glucose stock solution within 3 weeks.

9.3 Glucose-free Respiration Buffer:

Dissolve provided table in 50 mL of ddH₂O. Warm to assay temperature (usually 37°C), adjust to pH 7.4 with HCl and/or KOH and filter sterilize using a 0.22 µm filter. Store at 4°C. Use buffer within 3 weeks.

9.4 Glucose Oxidase Signal Control (GOx, 112.5 µg):

Reconstitute Glucose Oxidase in 40 µL of ddH₂O to generate a 3 mg/mL GOx stock solution. Aliquot into three equal volumes (10 µL) and store at -20°C. Each aliquot is sufficient for 2 wells (5 µL GOx/well).

9.5 Oligomycin (10.4 µg):

Reconstitute oligomycin in 95 µL of DMSO to generate a 100 µM (100X) stock solution. Aliquot into three equal volumes (30 µL) and store at -20°C. Each aliquot is sufficient for 30 wells (1 µM final concentration).

10. Sample Preparation

- Prepare a cell titration experiment to identify a suitable cell density for a specific cell type and conditions.
- Some cell lines (such as some glioblastoma cell lines) require the addition of glutamine to achieve their full glycolytic activity. This can be determined experimentally by adding Glutamine to the Respiration Buffer (typical concentration: 2 mM).
- 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 the acidification rate measured.
- If edge effects are observed, we recommend incubating the plate at RT for 30 minutes after plating before returning plate to cell culture. Alternatively, omit the outer wells of the plate.
- We recommend following the plate map suggested in figure 2 to facilitate assay procedure.

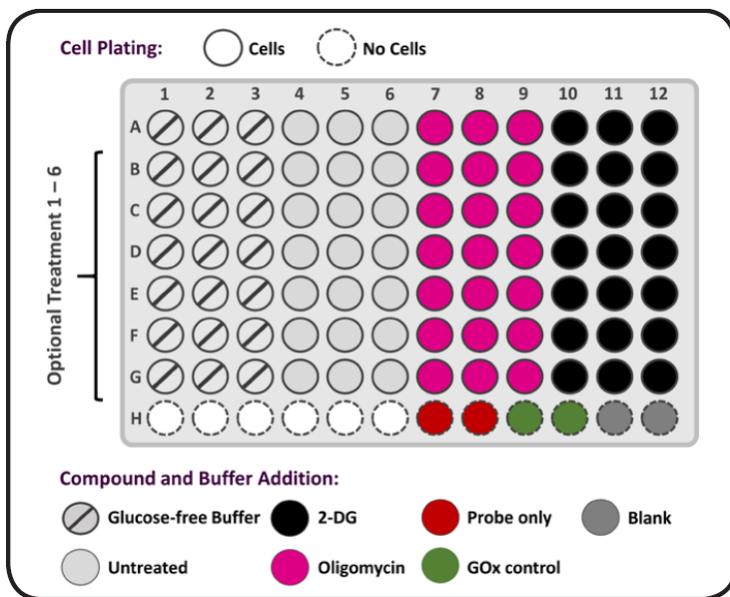


Figure 2. Recommended plate map for stress test assay with optional different treatments. Using this plate map, up to 7 different conditions can be tested simultaneously.

10.1 Cell plating:

10.1.1 Count cells and adjust to the desired plating density in culture medium – initial recommendation: $3-8 \times 10^4$ cells/well.

10.1.2 Return plate to incubator and culture overnight (~ 14 hours).

Δ Note: We recommend following plate map described in Figure 2 and leave row H with no cells.

Δ Note: If cells are cultured in CO₂ incubator overnight, you will need to place cells in a CO₂-free humidified incubator 2 hours prior to performing the assay (Step 11.2).

10.2 Compounds and assay controls set up:

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

- Oligomycin is typically used at 1 μM. However, the optimal concentration can be cell type dependent. Optimum concentrations can be determined by measuring ECA at varying oligomycin concentrations.
- Under standard measurement conditions, 50 mM 2-DG is usually sufficient to completely inhibit ECA caused by glycolysis. However, in some instances, higher concentrations may be necessary. Optimum concentrations can be determined by measuring ECA at varying 2-DG concentrations. Optional additional controls:
 - Glucose Oxidase is used as positive signal control for the assay.
- Compounds are typically added immediately prior to measurement. Longer times can be used as required: in this instances, compound should be present in both culture media (during incubation) and measurement media.
- If using test compounds, dilute so that the final concentration does not exceed 0.5% (v/v) and that volume added to well does not exceed 5 μL.

11. Assay Procedure

- This assay is designed to be assayed as a companion kit together in combination with Glycolysis Assay [Extracellular acidification] (ab197244). The Glycolysis 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 to running a Glycolysis Stress assay.
- We recommend that you assay all controls and samples in triplicate.
- Prepare all reagents 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). Ensure Respiration buffers are at the appropriate pH and pre-warmed at assay temperature (typically 37°C).

11.1 Prepare additional reagents:

- 11.1.1 Oligomycin 10X working stock (1 μ M end concentration): dilute 100X Oligomycin stock 1:10 in Respiration Buffer.
- 11.1.2 Glucose-containing Respiration Buffer: add Glucose Solution (Step 9.2) to Glucose-free Respiration Buffer to obtain desired glucose concentration (typically 5-20 mM). Buffer can be stored at 4°C for up to 3 weeks.

Δ Note: The user can define glucose concentration in the media to perform measurements at physiologically relevant glucose concentrations. Additionally, dependence of glucose concentration of glycolytic rate can be determined by measuring the ECA in the presence of oligomycin (ATP-limited glycolytic capacity) across a range of glucose concentrations.

- 11.1.3 Glycolysis Assay Reagent: prepare reagent as described in Glycolysis Assay [Extracellular acidification] (ab197244) user manual.

11.2 Wash cells:

Δ Note: Place cells in a CO₂-free humidified incubator 2 hours prior to performing the assay if cells were cultured in a CO₂ incubator overnight.

Δ Note: We recommend following plate map described in Figure 2 and leave row H with no cells.

11.2.1 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).

11.2.2 With a multichannel or repeater pipette, add 100 μL of the pre-warmed Glucose-Free Respiration Buffer to each well.

11.2.3 Repeat wash step one more time.

11.3 Add respiration buffer to wells:

- Glucose-free control (wells A-H 1-3) = 90 μL pre-warmed Glucose-free Respiration Buffer
- Untreated sample (wells A-H 4-6) = 90 μL pre-warmed Glucose-containing Respiration Buffer
- ATP-demand limited glycolysis (wells A-G 7-9) = 90 μL pre-warmed Glucose-containing Respiration Buffer
- Non-Glycolytic ECA (wells A-G 10-12) = 90 μL pre-warmed Glucose -containing Respiration Buffer
- Probe only (wells H7 and H8) = 90 μL pre-warmed Glucose -containing Respiration Buffer
- GOx Signal control (wells H9 and H10) = 90 μL pre-warmed Glucose -containing Respiration Buffer
- Blank control (wells H11 and H12) = 90 μL pre-warmed Glucose -containing Respiration Buffer

11.3.1 Add 10 μL of Glycolysis Assay Reagent (Step 11.1.2) to each well except to Blank Controls.

11.3.2 Add 10 μL of Glucose-containing Respiration Buffer to Blank Control wells (H11 and H12).

Δ Note: If measuring a full 96-well plate, we recommend diluting reconstituted ab197244 stock 1 in 10 in the relevant measurement media and, using a multichannel pipette, to add 100 μL to each well. Add 100 μL of Glucose-containing Respiration Buffer (no ab197244 reagent) to the Blank Control wells.

11.4 Treat cells:

Add relevant treatment or compounds to cells (see Section 12 for more detailed information on how to perform assays).

- 11.4.1 Add test compound or vehicle (typically 1-5 μL) to test wells (rows B-H, labeled "Optional Treatment 1-6" in Figure 2).
- 11.4.2 Add 10 μL oligomycin 10X Working Stock (Step 11.1.1) to wells to measure ATP-demand limited glycolysis (wells A-H 7-9).
- 11.4.3 Add 10 μL 2-DG Working Stock (Step 9.1) to wells to measure on-glycolytic ECA (wells A-H 10-12).
- 11.4.4 Add 10 μL Glucose-free Respiration Buffer to Glucose-free control (wells A-H 1-3) and Probe only (wells H7 and H8).
- 11.4.5 Add 5 μL GOx + 5 μL Glucose Solution (Step 9.2) to GOx Signal Control (wells H9 and H10).

11.5 Measurement:

- 11.5.1 Read plate immediately in a fluorescence microplate reader as described in the user manual instructions for Glycolysis Assay (ab197244).

11.6 Calculations:

- 11.6.1 Process data as described in Section 14 of the user manual instructions for Glycolysis Assay (ab197244).

12. Typical Assays/Data

12.1 Sample Glycolytic Characterization

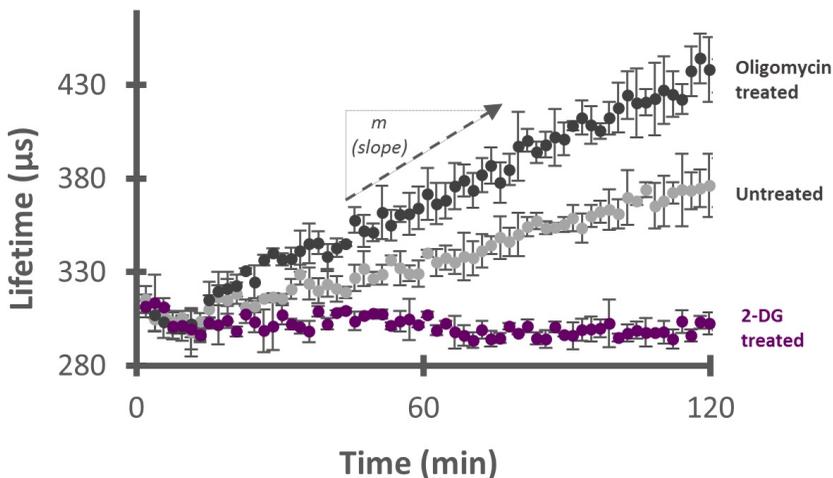


Figure 3. Extracellular acidification of A549 cells treated with 2-DG (purple) and oligomycin (black).

Basal Glycolysis: untreated cells (gray) show a steady increase in extracellular acidification caused by the export of glycolysis-derived protons. Basal glycolysis under given measurement conditions is dependent on factors such as glucose availability, demand for ATP and glycolytic intermediates, and the balance between oxidative phosphorylation and glycolytic ATP production.

Glycolytic Capacity: oligomycin treatment (black) inhibits F_1/F_0 ATPase, thereby preventing oxidative ATP generation and forcing the cells to compensate by increasing glycolytic flux to meet cellular ATP demand. The level of **Compensatory Glycolysis** can be determined by subtracting the Basal rate from the oligomycin-treated plate (Figure 4), This increase will be dependent on cellular ATP demand and the maximal glycolytic capacity under measurement conditions.

In instances where the ATP demand meets or exceeds the capacity of the cell to generate ATP glycolytically, the oligomycin-treated rate is reflective of *Maximal Glycolytic Capacity*. In instances where ATP demand does not exceed the capacity of the cell to generate ATP glycolytically, oligomycin-treated rate reflects *ATP-demand limited glycolysis* and informs of cellular response to the treatment condition.

Non-glycolytic ECA: 2-DG treatment (purple) inhibits glycolysis through competitive hexokinase inhibition allowing the measurement of non-glycolytic ECA; the acidification that is independent of transformation of glucose to lactate (eg, degradation of glycogen).

After conversion of the fluorescence lifetime signal into acidification rate (refer to Glycolysis Assay (ab197244) user manual for details), the calculated slopes (m , measure for acidification rates) can be used to determine the rate of non-glycolytic ECA (2-DG treated), basal glycolysis, ATP-demand limited glycolysis and the compensatory glycolysis (Figure 4).

Non-glycolytic ECA = m_{2-DG}

Basal Glycolysis = $m_{untreated} - m_{2-DG}$

ATP-demand limited glycolysis = $m_{Oligomycin}$

Compensatory Glycolysis = $m_{oligomycin} - m_{untreated}$

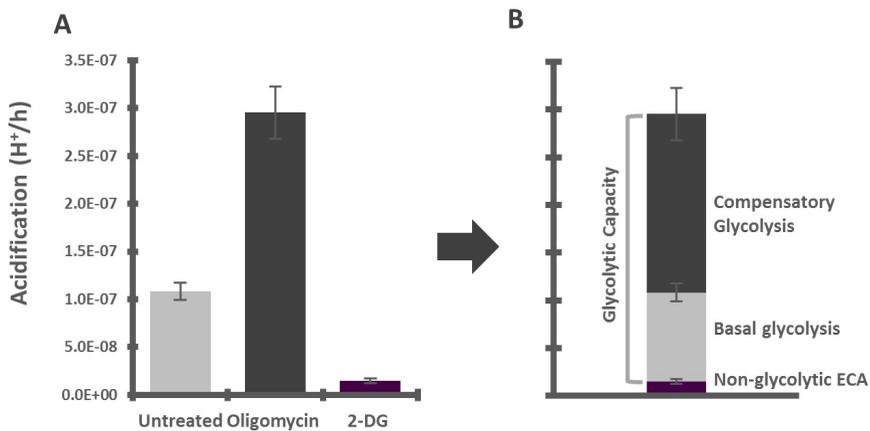


Figure 4. Glycolytic characterization of A549 cells.

12.2 Increasing ATP demand

As described in the previous section, the glycolytic rate observed upon oligomycin treatment is dependent on ATP demand and maximal glycolytic capacity under measurement conditions. To determine if oligomycin-treated glycolytic capacity is ATP-demand limited, cells can be pharmacologically challenged to increase ATP demand. In instances where increase ATP demand causes no further increase in measured ECA, this oligomycin-treated rate can be termed **Maximal Glycolytic Capacity**. Increased ECA in response to increased ATP demand illustrates that oligomycin-treated rates are ATP-limited.

Increased ATP demand can be induced by treatment with the ionophore monensin, which facilitates the electroneutral passage of monovalent cations across lipid membranes, thereby increasing the activity of the Na^+/K^+ -ATPase and by extension increasing ATP demand. Where cells have the capacity, compensatory glycolysis will increase to meet the increased ATP demand and additional detail on the glycolytic capacity of the cells under study can be obtained.

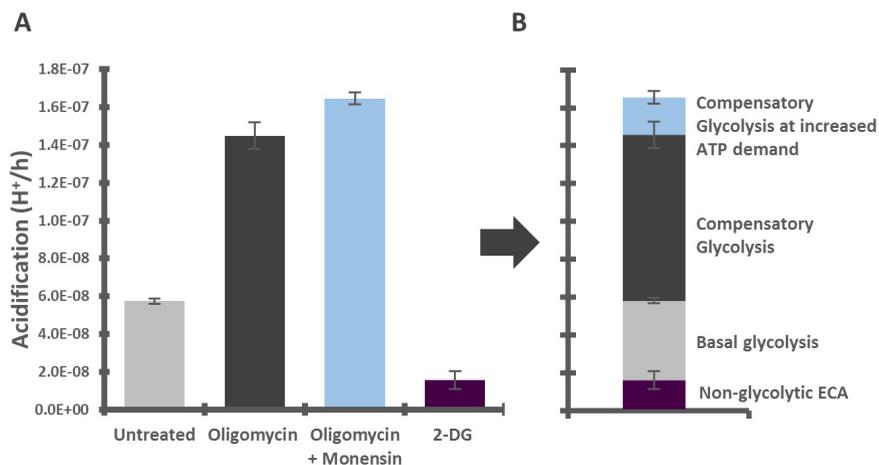


Figure 5. Increased compensatory glycolysis upon increased ATP demand in monensin-treated HepG2 cells.

13. FAQs

Q. Can I use other buffer than the Respiration Buffer provided in the kit?

A. The Respiration Buffer provided has been optimized for the assay. It contains 1 mM Potassium phosphate, 70 mM NaCl, 50 mM KCl, 0.8 mM MgSO₄, 2.4 mM CaCl₂.

Alternative media such as unbuffered DMEM can also be used for the assay, as long as care is taken to ensure a minimal buffering capacity.

14. Notes

Version 11 Last updated 6 January 2020

ab197244 Glycolysis Assay [Extracellular acidification]

For the measurement of extracellular acidification [ECA/ECAR] in cell populations and 3D culture models.

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

Table of Contents

15. Overview	25
16. Protocol Summary	26
17. Precautions	27
18. Storage and Stability	27
19. Limitations	28
20. Materials Supplied	28
21. Materials Required, Not Supplied	29
22. Technical Hints	30
23. Reagent Preparation	31
24. Plate Reader Set-Up	32
25. Signal Optimization	37
26. Assay Procedure	38
27. Assay Procedure for 384 well plate	40
28. Calculations	42
29. Typical data	43
30. Assay Throughput and Performance	44
31. Additional Assays/Data	45
32. FAQ	47
33. Notes	49

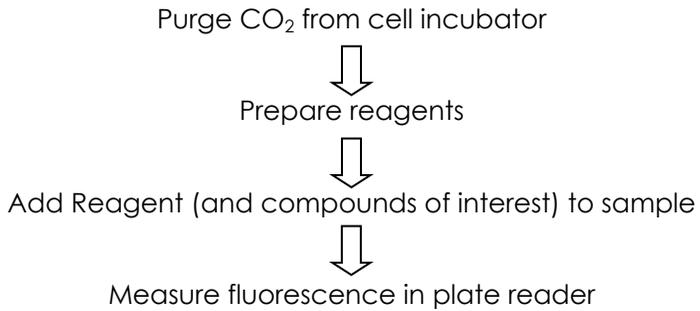
15. Overview

Glycolysis Assay [Extracellular Acidification] (ab197244) is an easy mix-and-measure, 96 or 384 well fluorescence plate reader-based approach for the analysis of extracellular acidification (ECA/ECAR). As lactate production is the main contributor to this acidification, Glycolysis Assay (ab197244) is a convenient and informative measuring tool of cellular glycolytic flux. Such measurements offer an important insight into the central role played by altered glycolytic activity in a wide array of physiological and pathophysiological processes, including cellular adaptation to hypoxia and ischemia, and the development and progression of tumorigenesis.

The pH-sensitive reagent is chemically stable and inert, water-soluble and cell impermeable. It exhibits a positive signal response (increased signal with increased acidification) across the biological range (pH 6 – 7.5). This performance, coupled with its spectral characteristics, make this kit the ideal choice for flexible, high-throughput assessment of extracellular acidification, overcoming many of the problems associated with the more cumbersome potentiometric pH approach. Rates of extracellular acidification are calculated from changes in fluorescence signal over time and, as the measurement is fully reversible, measurement of time courses and multiple drug treatments are possible.

The flexible plate reader format, allows multiparametric or multiplex combination with other products within this range. For example, this product in combination with the Extracellular O₂ Consumption Assay (ab197243) or the Extracellular O₂ probe (ab197242) allows the simultaneous real-time measurement of mitochondrial respiration and glycolysis and analysis of the metabolic phenotype of cells and the shift (flux) between the two pathways under pathological states.

16. Protocol Summary



17. 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 handled 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.

18. Storage and Stability

Store Reagent at 4°C and Buffer tablet at RT 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.

Δ Note: Reconstituted reagent is stable for 1 month.

19. 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.

20. Materials Supplied

Item	Quantity	Storage temperature (before prep)	Storage temperature (after prep)
Glycolysis Assay Reagent (lyophilized)	1 vial	4°C	-20°C
Respiration Buffer	1 tablet	RT*	4°C

*Store Respiration Buffer tablet at RT in dry/dark place (desiccant recommended).

21. 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 Table 1 in Instrument and Measurement Settings section for suitable plate readers
 - Double distilled water (ddH₂O)
 - 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
 - Sterilizing filter 0.22 μm
 - CO₂-free incubator
- For cells:
- Cell culture medium

22. 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 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 plates give optimal signal-to-noise ratios.
- For first time users, we recommend performing a Signal Optimization Step (see Section 11).

23. Reagent Preparation

Briefly centrifuge small vials at low speed prior to opening.

23.1 Respiration Buffer:

Reconstitute Respiration Buffer tablet in 50 mL of ddH₂O. Warm to assay temperature (usually 37°C), adjust to pH 7.4 and filter sterilize using a 0.22 µm filter. Store at 4°C in the dark.

23.2 Glycolysis Assay Reagent (lyophilized):

Reconstitute Assay Reagent in 1 mL of Respiration Buffer by gently aspirating and mixing 3-4 times.

Aliquot reagent so that you have enough volume to perform the desired number of assays. Store at -20°C in the dark.

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

24. Plate Reader Set-Up

24.1 Measurement Parameters

The Glycolysis Assay Reagent is a chemically stable and inert, water-soluble, cell impermeable pH-sensitive fluorophore.

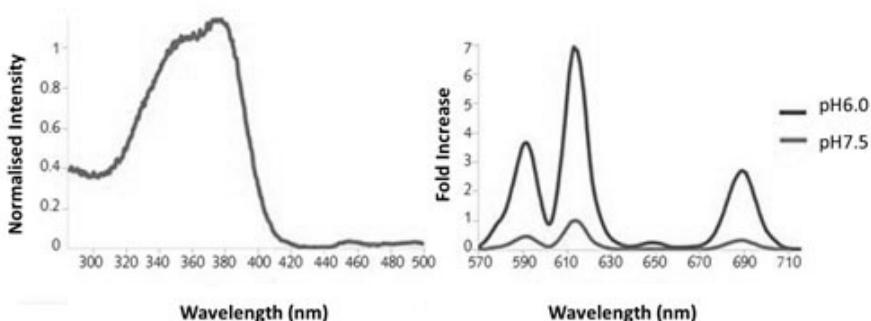


Figure 1. Excitation and emission spectra of the Glycolysis Assay Reagent. Left panel shows normalized excitation (Ex = 340-410 nm; Peak 360-380 nm). Right panel shows emission maxima fold increase (Em = 590, 615 and 690 nm) between pH 6.0 and pH 7.5.

- We strongly recommend using only fluorescence plate readers equipped with temperature control.
- This assay can also be used in non – TR-F Intensity mode on some plate readers, although we recommend running the Signal Optimization protocol (Section 12) to confirm an acceptable S:B, as well as optimizing cell seeding density (Figure 4).

24.2 Fluorescence measurements

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

24.2.1 Standard: TR-F Measurement

Time-resolved fluorescence (TR-F) measurement can be done in a wide range of commonly available plate readers (although users may see better performance using filter-based plate readers).

Optimal delay time is $\sim 100 \mu\text{s}$ and gate (integration) time is $100 \mu\text{s}$.

Δ Note: Glycolysis reagent should return Signal to Blank ratio (S:B) ≥ 3 .

24.2.2 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 and to express ECA as a function of $[\text{H}^+]$.

Δ Note: Glycolysis reagent should return Signal to Blank ratio (S:B) ≥ 10 .

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: $100 \mu\text{s}$ delay (D1), $30 \mu\text{s}$ measurement time (W1)
- Integration window 2: $300 \mu\text{s}$ delay (D2), $30 \mu\text{s}$ measurement time (W2)

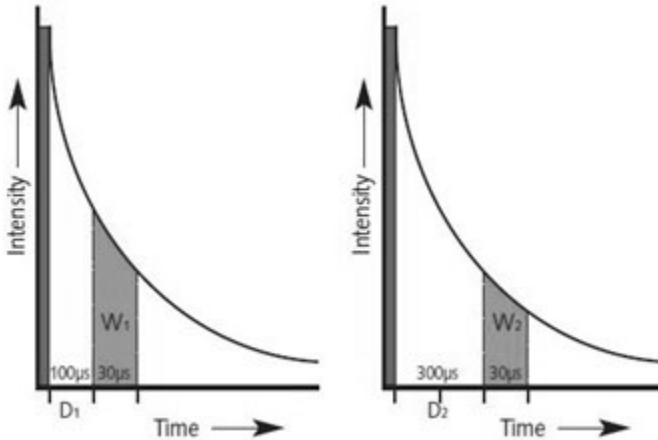


Figure 2. Illustrating dual read TR-F measurement.

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

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

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).

Δ Note: S:B for Integration window 2 is recommended to be ≥ 10 to allow accurate Lifetime calculation. Range of Lifetime values should be $\sim 200 \mu\text{s}$ for cells assayed in respiration buffer at approx. pH 7.4, increasing up to $> 400 \mu\text{s}$ upon acidification, and should only be calculated from samples containing reagent. Lifetime values should not be calculated from blank wells.

24.3 Instrument and Measurement Settings

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 μ s and 100 μ s.

Table 1. Recommended Instrument and Measurement Settings

Instrument	Optical Configuration	Intg1 (D ₁ /W ₁) Intg2 (D ₂ /W ₂)	Optimum Mode	Ex (nm) Em (nm)
BioTek: Synergy Neo, H4, 2 Cytation 3/5	Filter-based Top or bottom read	100 / 30 μ s 300 / 30 μ s	Dual-read TR-F (Lifetime)	Ex 360 \pm 40nm Em 620 \pm 10nm
BMG Labtech: CLARIOstar	Filter-based Top or bottom read	100 / 30 μ s 300 / 30 μ s	Dual-read TR-F (Lifetime)	Ex 340 \pm 50nm (TR-EX L) Em 615 \pm 10nm (BP-615)
BMG Labtech: FLUOstar Omega / POLARstar Omega	Filter-based Top or bottom read	100 / 30 μ s 300 / 30 μ s	Dual-read TR-F (Lifetime)	Ex 340 \pm 50nm (TR-EXL) Em 615 \pm 10nm (BP-615)
BMG Labtech: PHERAstar FS	Filter-based Top read (HTRF Module)	100 / 30 μ s 300 / 30 μ s	Dual-read TR-F (Lifetime)	Ex 337 nm Em 620 nm
Perkin Elmer: VICTOR series/ X4, X5	Filter-based Top read	100 / 30 μ s 300 / 30 μ s	Dual-read TR-F (Lifetime)	Ex 340 \pm 40nm (D340) Em 615 \pm 8.5nm (D615)

Instrument	Optical Configuration	Intg1 (D ₁ /W ₁) Intg2 (D ₂ /W ₂)	Optimum Mode	Ex (nm) Em (nm)
Perkin Elmer: EnVision	Filter-based Top read	100 / 50µs 300 / 50µs	Dual-read TR-F (Lifetime)	Ex 340 ±60nm (X340) Em 615 ± 8.5nm (M615)
Tecan: Infinite M1000 Pro/ F200 Pro	Monochromator Filter-based Top or bottom read	100 / 100µs 300 / 30 µs	Dual-read TR-F (Lifetime)	Ex 380±20nm Em 615±10nm
BMG Labtech: FLUOstar Optima / POLARstar Optima	Filter-based Top or bottom read	100 / 100µs n/a	TR-F	Ex 340 ± 50nm (TR-EXL) Em 615 ± 10nm (BP-615)
Perkin Elmer: EnSpire	Monochromator Top or bottom read	100 / 100µs n/a	TR-F	Ex 380 nm Em 615 nm
Tecan: Infinite M200 Pro/ Sapphire / Genios Pro	Monochromator Filter-based / Top or bottom read	100 / 100µs n/a	TR-F	Ex 380±20nm Em 615±10nm
Mol. Devices: SpectraMax / Flexstation / Gemini	Monochromator based Top or bottom read	50 /250µs n/a	TR-F	Ex 380nm Em 615nm

25.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.

25.1 Signal Optimization:

- 25.1.1 Prepare 8 replicate wells of a 96-well plate, by adding 150 μ L pre-warmed culture medium to each well (A1-A4, B1-B4).
- 25.1.2 Add 10 μ L reconstituted Glycolysis Assay Reagent to 4 of the replicate wells (A1-A4) and 10 μ L ddH₂O, PBS or media to the remaining replicates wells (B1-B4).
- 25.1.3 Read plate immediately in a fluorescence plate reader over 30 minutes (read every 2-3 minutes).
- 25.1.4 Examine Signal Control well (A1-A4) and Blank Control well (B1-B4) readings (linear phase) and calculate Signal to Blank (S:B) ratio.

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

	1	2	3	4
A	Respiration Buffer + Glycolysis Reagent			
B	Respiration Buffer	Respiration Buffer	Respiration Buffer	Respiration Buffer

26. 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).
- The amount of signal change will be directly dependent on the rate of glycolytic flux of the cell type being measured. We recommend using a medium – to – high cell density per well as a starting point, and reducing cell numbers as required.
- 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
Glucose Oxidase (positive signal control)	1 mg/mL in ddH ₂ O
FCCP (ETC uncoupler) (positive control, increases ECA)	Titration recommended to establish best concentration
Oxamic acid (negative control, decreases ECA)	750 mM in ddH ₂ O

26.1 Pre-assay preparation – Adherent cells:

26.1.1 Seed cells in a 96-well plate at a density of $3 - 8 \times 10^4$ cells/well in 200 μ L culture medium.

26.1.2 Incubate overnight in a CO₂ incubator at 37°C.

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

Where cells are cultured in CO₂ incubator overnight, it is important to purge the media and plasticware of CO₂ prior to conducting the assay as residual CO₂ may contribute to acidification. Perform a CO₂ purge by incubating cells in a CO₂-free incubator at 37°C with 95% humidity, approx. 3 hours prior to performing the Glycolysis assay measurement.

26.2 Pre-assay preparation – Suspension cells:

- 26.2.1 Harvest suspension cells.
- 26.2.2 Wash cells in Respiration Buffer.
- 26.2.3 Seed cells in a 96-well plate at a density of $2.5 - 5 \times 10^5$ cells/well in 150 μ L Respiration Buffer.
- 26.2.4 Proceed to plate relevant controls (Step 12.3.4).

26.3 Plate loading:

- 26.3.1 Remove spent culture media from all assay wells and wash cells with 100 μ L of Respiration Buffer.
- 26.3.2 Repeat washing step.
- 26.3.3 Add 150 μ L of Respiration Buffer to all wells containing cells.
- 26.3.4 Blank controls (we suggest using wells H11 and H12): add 150 μ L of Respiration Buffer.
- 26.3.5 Optional positive control: add 10 μ L of glucose oxidase solution (1 mg/mL in ddH₂O) to wells containing 150 μ L Respiration Buffer only.
- 26.3.6 Optional negative control: add 10 μ L of oxamic acid solution (750 mM in ddH₂O) to wells containing cells.

26.4 Assay set up:

- 26.4.1 Add 10 μ L reconstituted Glycolysis Assay Reagent to each sample well and positive/negative control wells.
- 26.4.2 Add 10 μ L of Respiration Buffer to blank control wells.
- 26.4.3 Add 1 – 10 μ L test compound (vehicle control and/or stock) to the wells.

Δ Note: We recommend keeping the volume of added compound low to minimize any potential effects of solvent vehicle.

Δ Note: Plate preparation time should be kept to a minimum.

26.5 Measurement:

- 26.5.1 Insert the prepared plate into a fluorescence plate reader pre-set to the measurement temperature (typically 37°C).
- 26.5.2 Measure Glycolysis assay signal at 1.5 min intervals for ≥ 120 minutes using excitation and emission wavelengths of Ex/Em = 380/615 nm respectively (see Table 1 for Instrument and Measurement Setting table for instrument specific information).

27. 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 12.

27.1 Pre-assay preparation – Adherent cells:

27.1.1 Seed cells in a 384-wp at a density of $2 - 4 \times 10^4$ cells/well in 100 μL culture medium.

27.1.2 Incubate overnight in a CO_2 incubator at 37°C .

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

Δ Note: Where cells are cultured in CO_2 incubator overnight, it is important to purge the media and plasticware of CO_2 prior to conducting the assay as residual CO_2 may contribute to acidification. Perform a CO_2 purge by incubating cells in a CO_2 -free incubator at 37°C with 95% humidity, approx. 3 hours prior to performing the Glycolysis assay measurement.

27.2 Pre-assay preparation – Suspension cells:

27.2.1 Harvest suspension cells.

27.2.2 Wash cells in Respiration Buffer.

27.2.3 Seed cells in a 384-wp at a density of $1.2 - 2.5 \times 10^5$ cells/well in 75 μL Respiration Buffer.

27.2.4 Proceed to plate relevant controls (Step 13.3.4).

27.3 Plate loading:

27.3.1 Remove spent culture media from all assay wells and wash cells with 50 μL of Respiration Buffer.

27.3.2 Repeat washing step.

27.3.3 Add 75 μL of Respiration Buffer to all wells containing cells.

27.3.4 Blank controls (we suggest using wells H11 and H12): add 75 μL of Respiration Buffer.

27.3.5 Optional positive control: add 5 μL of glucose oxidase solution (1 mg/mL in ddH₂O) to wells containing 75 μL Respiration Buffer only.

27.3.6 Optional negative control: add 5 μL of oxamic acid solution (750 mM in ddH₂O) to wells containing cells.

27.4 Assay set up:

- 27.4.1 Add 5 μ L reconstituted Glycolysis Assay Reagent to each sample well and positive/negative control wells.
- 27.4.2 Add 5 μ L of Respiration Buffer to blank control wells.
- 27.4.3 Add 1 – 5 μ L test compound (vehicle control and/or stock) to the wells.

Δ Note: we recommend keeping the volume of added compound low to minimize any potential effects of solvent vehicle.

Δ Note: plate preparation time should be kept to a minimum.

27.5 Measurement:

- 27.5.1 Insert the prepared plate into a fluorescence plate reader pre-set to the measurement temperature (typically 37°C).

Measure Glycolysis assay signal at 1.5 min intervals for ≥ 120 minutes using excitation and emission wavelengths of Ex/Em = 380/615 nm respectively (see Table 1 for Instrument and Measurement Setting table for instrument specific information).

28. Calculations

- Plot the Blank control well-corrected Glycolysis Assay Intensity or Lifetime values versus Time (min).
- Select the linear proportion of the signal profile (avoiding any initial lag or subsequent plateau) and apply linear regression to determine the slope (ECA) and correlation coefficient for each well.

Δ Note: this approach is preferable to calculating a slope from averaged profiles.

- 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.

29. Typical data

Data provided for **demonstration purposes** only.

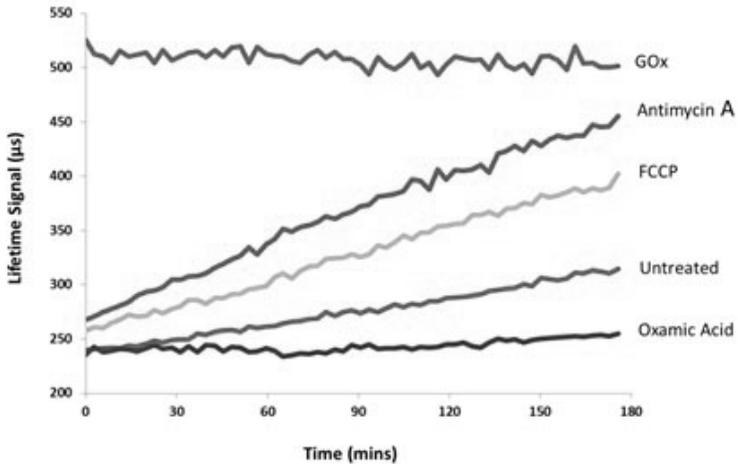


Figure 3. Typical Lifetime profile of Glycolysis Assay for adherent cells (HepG2), treated with modulator compounds, including oxamic acid (recommended negative control as inhibits conversion of pyruvate to lactate). Effect of glucose oxidase (GOx) as positive signal control is illustrated schematically.

30. Assay Throughput and Performance

Figure 4A illustrates a typical read out of showing the parallel analysis of 96 individual samples. A serial dilution of HepG2 is presented in Figure 4B, with increasing numbers cell numbers causing an increased rate of acidification. This is seen as an increased rate of signal change.

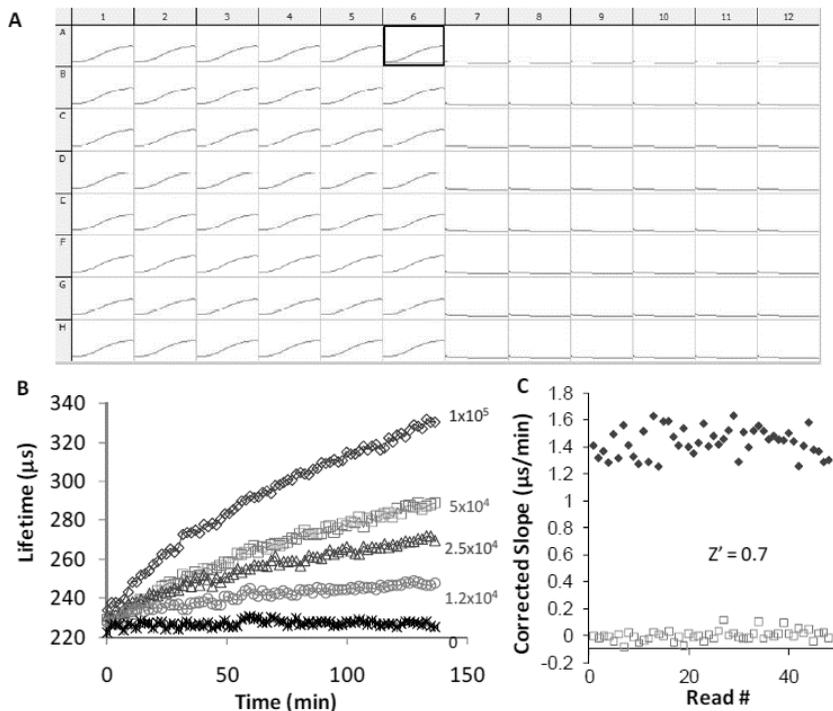


Figure 4. Analysis of cellular respiration of HepG2 cells. (A) 96-wp based analysis of extracellular acidification (cell profiles on left, control proliferates on right). (B) Acidification profiles for HepG2 cultured in L15 medium at the indicated seeding concentration. (C) Z' factor analysis for HepG2 plated in L15 medium at $\sim 10^5$ cells/well. Data generated on FLUOstar Omega (A) and on a Victor X4, PerkinElmer (B & C).

The performance of the assay is highlighted in Figure 4C, with a coefficient of variance (%CV) below 5%. Z' factor analysis assesses assay performance in terms of signal window and measurement reproducibility and shows excellent performance ($Z' = \sim 0.7$).

31. Additional Assays/Data

31.1 Titration of cell density

To determine an optimal cell seeding density for performing this assay with new cell types, seed cells (in replicate) with a range of seeding densities (typically 0, 10^4 , 2×10^4 , 4×10^4 , 6×10^4 , 8×10^4 cells/well).

Plot the data generated as a function of intensity or Lifetime values versus time, as illustrated in Figure 4.

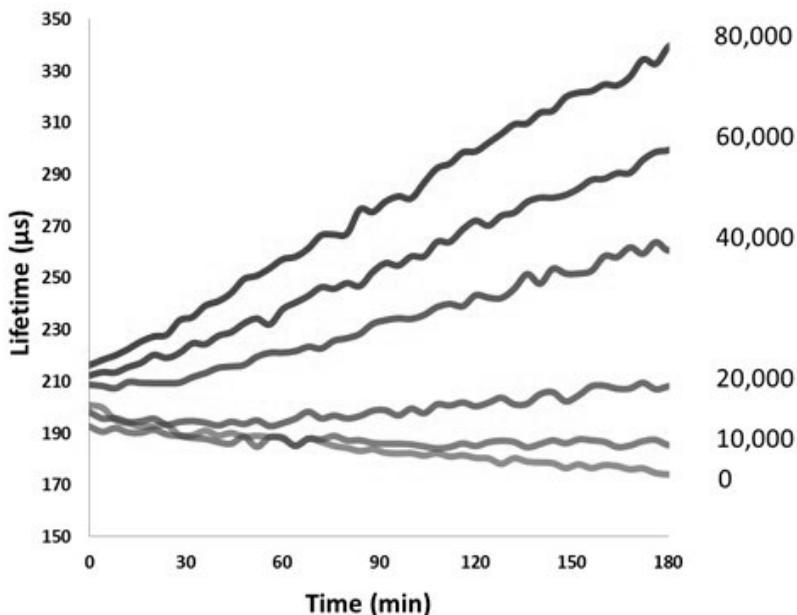


Figure 5. Extracellular acidification rate (ECA) profiles are shown for A549 cells seeded at 0, 10^4 , 2×10^4 , 4×10^4 , 6×10^4 , 8×10^4 cells/well. In this experimental example, a seeding density of 40,000 cells/well was chosen for further study as it provided a suitable balance between ECA response and cell availability.

31.2 Cellular energy flux analysis

Multiparametric (or multiplex) combination of Glycolysis Assay together with Extracellular O₂ Consumption Assay (ab197243) 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.

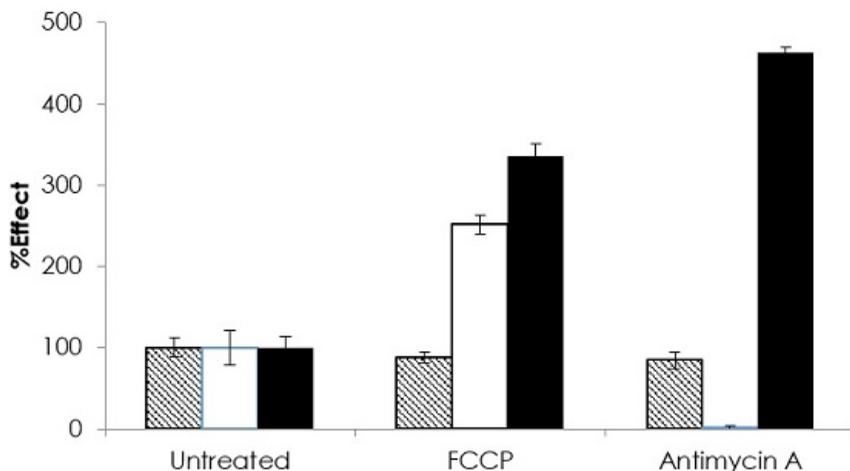


Figure 6. Cellular Energy Flux for HepG2 cells (seeded at 65,000 per well), treated with a combination of drug compounds modulating the ETC (Antimycin A [1 μ M] and FCCP [2.5 μ M]), shown as a percentage relative to untreated control cells. Comparative measurements were taken with Extracellular Oxygen 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)).

32. FAQ

Q. What can I do if Signal to Blank (S:B) ratio is not as high as expected?

A. Glycolysis Reagent should return a S:B ratio ≥ 3 .

The following options may be helpful to improve S:B if the ratio is not as high as expected:

- Increase Gain (PMT) setting or flash energy
- Adjust TR-F focal height
- Increase length of integration time, the same for both delay windows.
- Repeat as top or bottom-read, respectively.
- Increase volume of Glycolysis Reagent (15 μ l)
- Contact Instrument Supplier for further options

Q. What do I do if I cannot detect any signal in wells containing cells and Glycolysis Reagent (or I can detect a signal but the slope rate appears very low)?

A. There are few things you should check:

- Check you have the correct instrument settings.
- Perform Signal Optimization.
- Include GOx control (max signal).
- Increase cell density.
- Check pH of pre-warmed Respiration Buffer and correct as necessary, as pH can drift over time.

Q. What do I do if I can detect a signal in wells containing cells and Glycolysis Reagent, but the slope (rate) falls initially or is variable from well to well?

A. There are few things you should check:

- Check cell seeding and pipetting consistency.
- Increase cell density.
- Ensure plate, instrument and all culture media and stock solutions are pre-warmed at 37°C prior to use.
- Reduce plate preparation times.

Q. Can I use other buffer than the Respiration Buffer provided in the kit?

A. The Respiration Buffer provided in the kit has been optimized for the assay. It contains 1 mM Potassium Phosphate, 20 mM Glucose, 70 mM NaCl, 50 mM KCl, 0.8 mM MgSO₄, 2.4 mM CaCl₂.

Alternative media and supplements may be used as required (such as unbuffered DMEM), as long as care is taken to ensure a minimal buffering capacity.

33. Notes

Technical Support

Copyright © 2014-2020 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