

Glycolysis/OXPHOS Assay Kit

Technical Manual

General Information

Adenosine triphosphate (ATP), synthesized by glycolysis and mitochondrial oxidative phosphorylation, is a crucial energy source for living cells. Many cancer cells produce ATP by glycolysis. However, recent studies have suggested that if glycolysis is suppressed in cancer cells, their ATP synthesis shifts to oxidative phosphorylation by enhancing mitochondrial function^{1,2}. It is expected that understanding these phenomena will be helpful for elucidation of the mechanism of anticancer drugs and will lead to the development of therapeutic strategies for cancer, neurodegenerative disease, and various diseases.

The Glycolysis/OXPHOS Assay Kit enables analysis of cellular glycolytic capacity, whether cells rely on glycolysis or oxidative phosphorylation for energy production, and metabolism shift by using inhibitors (Oligomycin and/or 2-DG). The kit includes all the reagents required for these evaluations.

Kit Contents

Dye Mixture (Blue cap)	×1	Use for Lactate Assay
Lactate Standard (10 mmol/l) (Blue cap)	150 µl×1	
LDH Solution (Purple cap)	12 µl×1	
Lactate Assay Buffer	5.5 ml×1	
Reconstitution Buffer (White cap)	550 µl×1	
Substrate (Green cap)	×1	Use for ATP Assay
Luciferase Solution (Green cap)	10 µl×1	
ATP Assay Buffer	5.5 ml×1	
Oligomycin (Yellow cap)	×1	Use as an inhibitor
2-DG (Red cap)	140 µl×1	

※ This kit includes the reagents corresponding to 48 wells of a 96-well microplate. The number of samples that can be evaluated with this kit is posted on the product web page.

※ Evaluation with this kit is performed by relative evaluation using absorbance and/or luminescence values.

Storage Conditions

Store at 0–5 °C

Required Equipment and Materials

- Microplate reader (450 nm filter and luminescence)
- 96-well microplate, 96-well white microplate
- Incubator (37 °C)
- 20–200 µl multichannel pipette
- 100–1000 µl, 20–200 µl, 2–20 µl micropipettes
- Dimethyl sulfoxide (DMSO)
- Conical tube

Precautions

- Equilibrate reagents to room temperature prior to use.
- Briefly centrifuge the tubes containing Lactate Standard (blue cap), LDH Solution (purple cap), Reconstitution Buffer (white cap), Luciferase Solution (green cap) and 2-DG (red cap) before opening to remove all content from the walls of the tube and inside the cap.
- Pipette the LDH Solution (used for Lactate Assay) before use to ensure the enzyme is homogenous.
- Analysis of samples in triplicate is recommended for accuracy.
- Since the enzymatic reaction starts immediately after adding Lactate working solution to a well, use a multichannel pipette to minimize the experimental error from time lag in pipetting.
- Dye Mixture and Substrate are stored in a glass bottle with an aluminum cap. Handle with caution, wear gloves.

Preparation of Solution (inhibitor)

Preparation of Oligomycin stock solution (10 mmol/l)

Add 10 µl DMSO to the Oligomycin tube (yellow cap) and dissolve by pipetting.

※ It is difficult to see the Oligomycin reagent. After adding DMSO, repeat pipetting several times and mix well.

※ The Oligomycin stock solution is stable for 2 months when stored -20 °C.

Preparation of Solutions and Measurement (Lactate Assay)

Preparation of Dye Mixture stock solution

Add all Reconstitution Buffer (white cap) to a Dye Mixture vial. Close the cap and dissolve the contents completely.

※ Transfer the Dye Mixture stock solution to the vial of the Reconstitution Buffer and store it at 0–5 °C with protection from light.

Dye Mixture stock solution is stable for 4 months under these conditions.

Preparation of Lactate working solution

(1) Add Dye Mixture stock solution to a conical tube and dilute it with Lactate Assay Buffer.

(2) Add LDH Solution (purple cap) to the solution prepared in step (1).

※ Refer to Table 1.

※ Lactate working solution is light sensitive. Prepare the solution just before use and protect it from light by covering it with aluminum foil. Please use up Lactate working solution within that day.

Table 1. Examples of Lactate working solution preparation

	for 24 well	for 48 well
Dye Mixture stock solution	250 µl	500 µl
Lactate Assay Buffer	2.25 ml	4.5 ml
LDH Solution (Purple cap)	5 µl	10 µl

Sample preparation

Prepare cell culture supernatant sample.

※ Please prepare samples diluted with double-deionized H₂O (ddH₂O) so that the absorbance is lower than that of 1 mmol/l Lactate standard solution (see Figure 1).

※ In case a medium contains serum, read the blank absorbance (serum containing medium) as background control and subtract its value from the absorbance of each sample.

※ Required sample amount is 20 µl for each well.

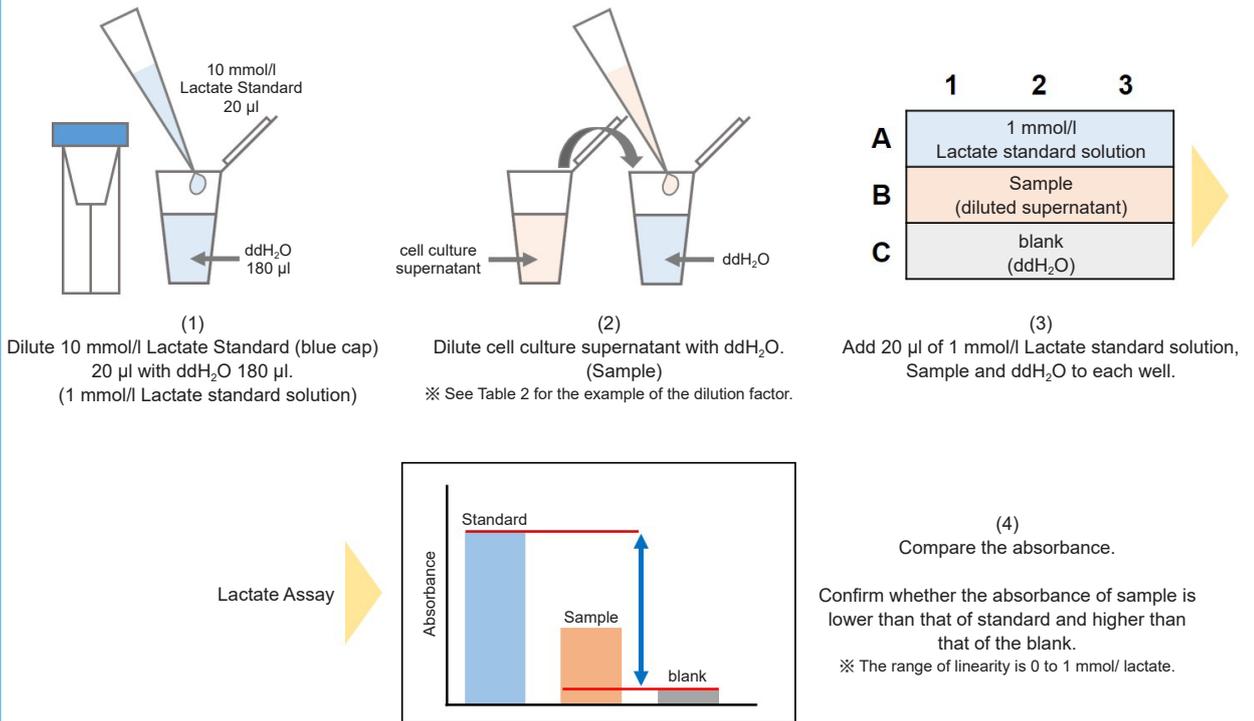
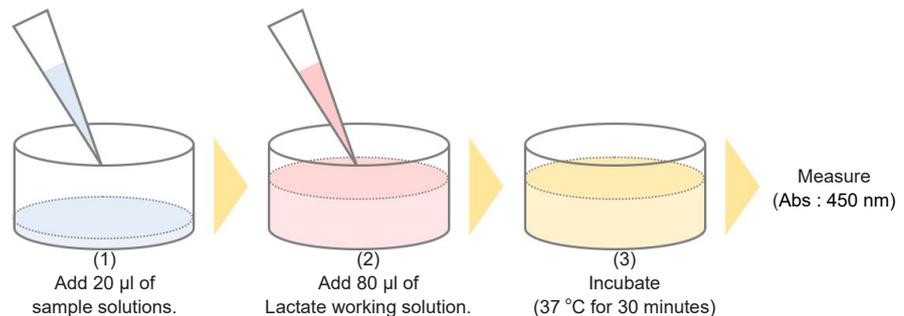


Figure 1. Method for dilution of samples for Lactate Assay

Table 2. Example of dilution

		Inhibitor treatment time		
		3 hours	5 hours	24 hours
24-well	HeLa, HepG2, A549 (5×10^4 cells/well)	10-fold	10-fold	20-fold
	WI-38 (5×10^4 cells/well)	10-fold	10-fold	15-fold
	Jurkat (1×10^5 cells/well)	10-fold	10-fold	15-fold
96-well	HeLa, HepG2, A549 (5×10^3 cells/well)		10-fold	15-fold
	WI-38 (5×10^3 cells/well)		10-fold	10-fold
	Jurkat (1×10^4 cells/well)		10-fold	10-fold

Measurement



(1) Add 20 µl of sample solutions to each well.

※ To obtain accurate data, triplicate measurements are recommended.

(2) Add 80 µl of Lactate working solution to each well.

※ Because the enzymatic reaction starts immediately after the Lactate working solution is added, use a multichannel pipette to minimize time lags in pipetting.

(3) Incubate the microplate at 37 °C for 30 minutes.

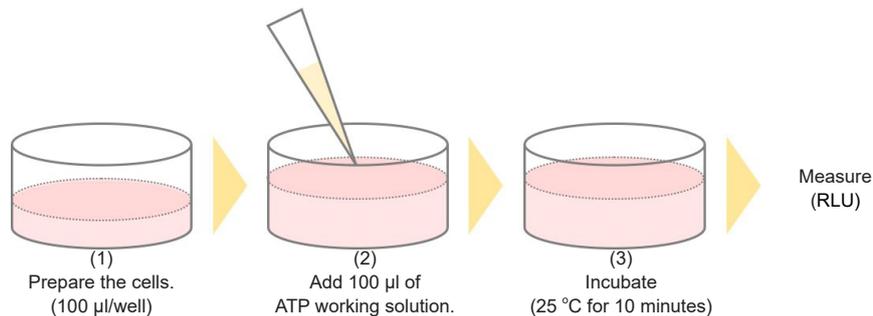
※ Seal the microplate during incubation to prevent evaporation.

(4) Measure the absorbance at 450 nm with a microplate reader.

Preparation of ATP working solution

- (1) Add 1 ml of ATP Assay Buffer to the Substrate vial.
 ※ Substrate vial is capped under vacuum pressure. Please open it carefully. Do not open quickly to avoid spilling contents.
- (2) Close the cap of the Substrate and dissolve the contents completely and transfer it to the ATP Assay Buffer bottle.
- (3) Add 10 µl of Luciferase Solution (green cap) to the solution prepared in step (2).
- (4) Mix by gently inverting the contents to obtain ATP working solution.
 ※ The ATP working solution is stable for 1 month when stored at -20 °C.

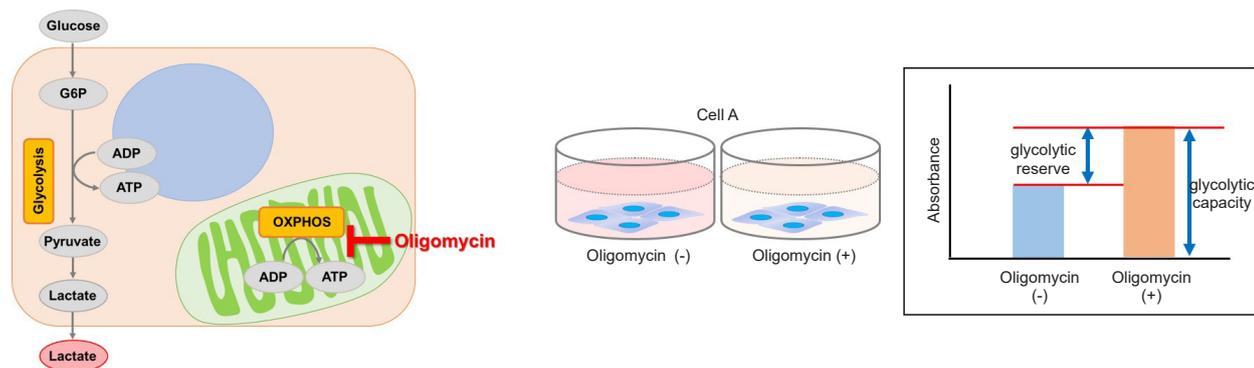
Measurement



- (1) Prepare the cells in 96-well white microplate.
 ※ To obtain accurate data, triplicate measurements are recommended.
- (2) Add 100 µl of ATP working solution to each well.
 ※ Mix the contents by using the process of reverse pipetting to avoid air bubbles that could influence the luminescent signal.
 ※ After adding the ATP working solution, mix the contents for 2 minutes on an orbital shaker or similar.
 ※ Please protect the plate from light by covering it with aluminum foil because the luminescent signal is light sensitive.
- (3) Incubate the microplate at 25 °C for 10 minutes in a microplate reader.
 ※ Please incubate the microplate in an incubator which has been temperature-settled at 25 °C or at a room temperature around 25 °C.
 ※ The above incubation is necessary to efficiently extract ATP from cells and for a stable luminescent signal.
- (4) Measure luminescence with a microplate reader.

Overview of Analysis

The purpose of this evaluation is to know the enhanced level of glycolysis in the cells using lactate production. To confirm the cellular glycolytic capacity, inhibit ATP synthesis by oxidative phosphorylation with a treatment using Oligomycin.



Reagents

Table 3. Reagents for analysis of cellular glycolytic capacity

Dye Mixture (Blue cap)	×1	Use for Lactate Assay
Lactate Standard (10 mmol/l) (Blue cap)	150 µl×1	
LDH Solution (Purple cap)	12 µl×1	
Lactate Assay Buffer	5.5 ml×1	
Reconstitution Buffer (White cap)	550 µl×1	
Oligomycin (Yellow cap)	×1	Use as an inhibitor

Preparation of Oligomycin solution

(1) Dilute the 10 mmol/l Oligomycin stock solution 1,000-fold with culture medium (to give 10 μ mol/l Oligomycin solution).

※ For example, to prepare 1 ml of Oligomycin solution, mix 1 μ l of 10 mmol/l Oligomycin stock solution and 1 ml of culture medium.

(2) Dilute the solution prepared in step (1) 8-fold with culture medium to give 1.25 μ mol/l Oligomycin solution.

※ Where cells are to be stimulated with a drug, add the drug to the medium used to prepare the Oligomycin solution so that the drug concentration is the same as required for stimulation (see Figure 2).

※ Use Oligomycin solution within the same day (it cannot be stored).

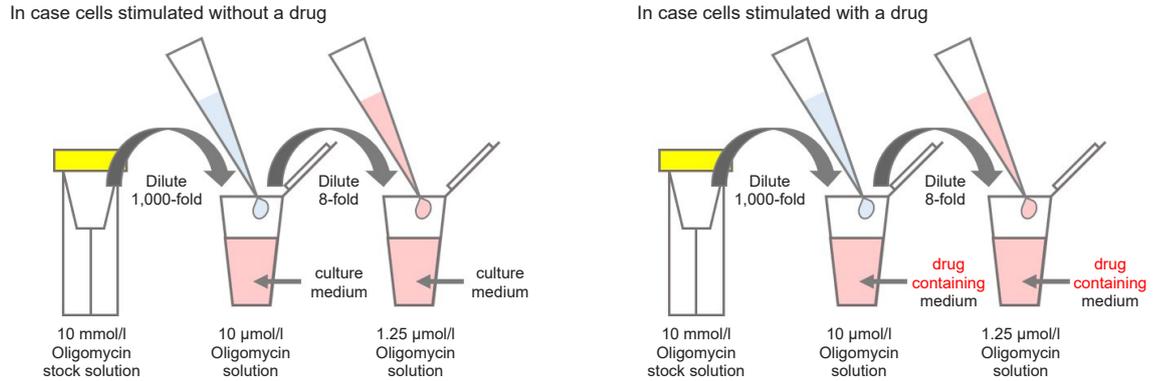
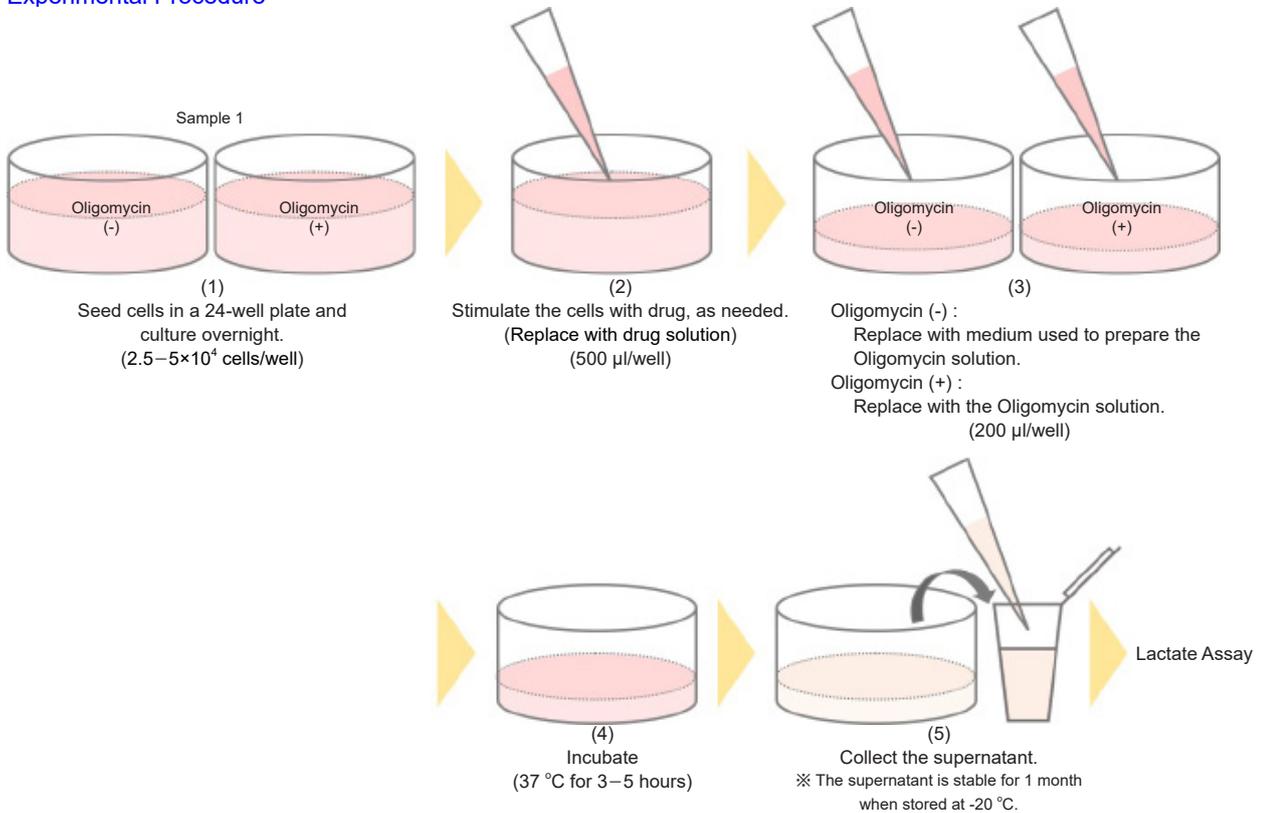


Figure 2. Preparation of Oligomycin-containing medium

Experimental Procedure



(1) Seed cells in a 24-well plate (2.5–5×10⁴ cells/well) and culture overnight in an incubator (37 °C, 5% CO₂).

※ Wells with and without Oligomycin treatment are required for each sample.

(2) Remove the medium and stimulate the cells with the drug, as needed.

※ Add drug solution to wells.

(3) Remove the medium and add 200 μ l of 1.25 μ mol/l Oligomycin solution to test sample Oligomycin (+) wells.

(4) Add 200 μ l of the medium used to prepare the Oligomycin solution to test sample Oligomycin (-) wells.

(5) Incubate the microplate for 3-5 hours in an incubator (37 °C, 5% CO₂).

(6) Collect 100 μ l of each cell culture supernatant into a 1.5 ml microtube and perform the Lactate Assay.

※ See Figure 3 for an assay plate format when performing the Lactate Assay.

(7) Compare the absorbance values of with and without Oligomycin treatment.

	1	2	3
A	Sample 1 Oligomycin (-)		
B	Sample 1 Oligomycin (+)		
C	Sample 2 Oligomycin (-)		
D	Sample 2 Oligomycin (+)		
E	medium only (background)		

Figure 3.

Example of an assay plate format (n=3)

Experimental Example : Change in glycolytic capacity of glycolytic inhibitor-treated HeLa cells

- (1) HeLa cells were seeded in four wells of a 24-well plate (2.5×10^4 cells/well in MEM containing 10% fetal bovine serum and 1% penicillin-streptomycin) and cultured overnight in an incubator (37 °C, 5% CO₂).
- ※ Two wells were prepared for glycolytic inhibitor stimulation and two wells for the control treatment.
- (2) The medium was removed, and glycolytic inhibitor solution (500 µl, in MEM containing 10% FBS) was added to two cell-seeded wells of the 24-well plate, and MEM containing 10% FBS (500 µl) was added to the other two cell-seeded wells.
- (3) The cells were cultured overnight in an incubator (37 °C, 5% CO₂).
- (4) The medium was removed, and 1.25 µmol/l Oligomycin solution (200 µl) was added to the Oligomycin (+) wells.
- (5) The medium (200 µl) used to prepare the Oligomycin solution was added to the Oligomycin (-) wells.
- (6) The cells were cultured for 3 hours in an incubator (37 °C, 5% CO₂).
- (7) After incubation, 100 µl of each cell culture supernatant was transferred to a 1.5 ml microtube and diluted 10-fold with ddH₂O to prepare sample solutions.
- ※ For background control, MEM only was also diluted 10-fold with ddH₂O.
- (8) The sample solutions (20 µl each) were added to individual wells of a 96-well microplate.
- (9) Lactate working solution (80 µl) was added to each well.
- (10) The assay plate was incubated at 37 °C for 30 minutes.
- (11) The absorbance at 450 nm was measured using a microplate reader, and the absorbance values of with and without Oligomycin treatments were compared.

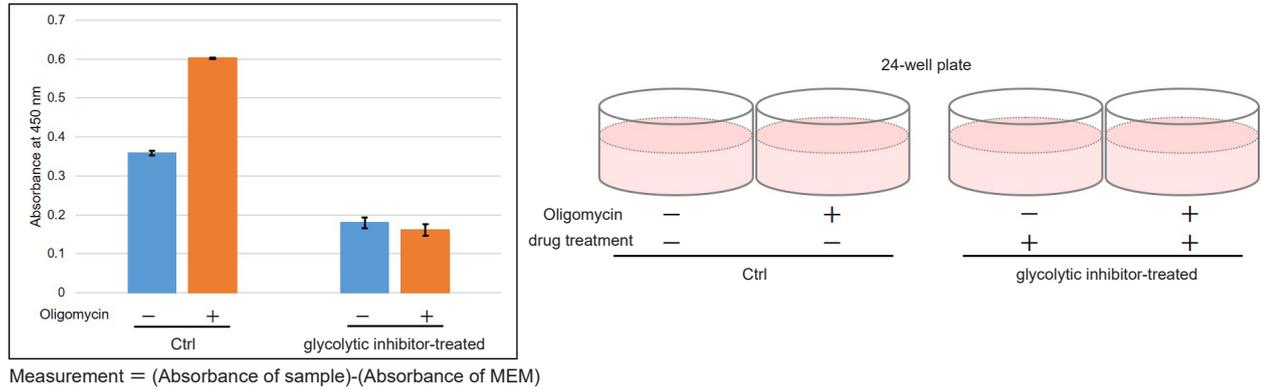
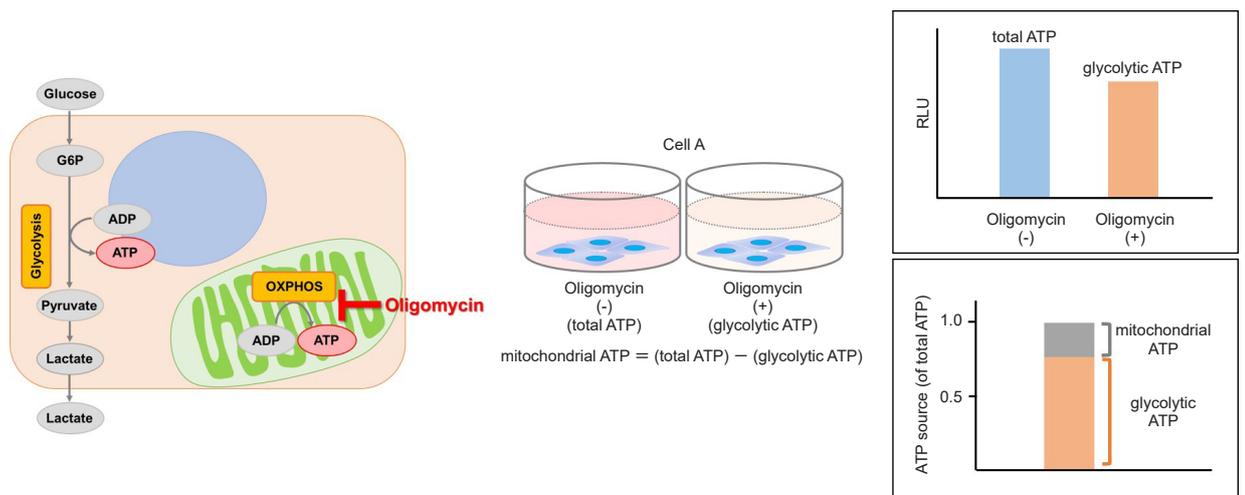


Figure 4. Comparison of glycolytic capacity of glycolytic inhibitor-treated and non-treated (Ctrl) HeLa cells
 In case HeLa cells are treated with Oligomycin to inhibit ATP synthesis by oxidative phosphorylation, glycolytic inhibitor non-treated (Ctrl) cells can enhance glycolysis, but glycolytic inhibitor-treated cells cannot enhance glycolysis (decreased glycolytic capacity).

General Protocol ②
(Analysis of Metabolism Shift)

Overview of Analysis

The values can be obtained with this evaluation allows you to know whether cells show a metabolic shift. The amounts of ATP produced by glycolysis and mitochondrial oxidative phosphorylation are determined from the amounts of ATP when the cells were treated with and without Oligomycin (which inhibits ATP synthesis in oxidative phosphorylation).



Reagents

Table 4. Reagents for analysis of metabolism shift

Substrate (Green cap)	×1	Use for ATP Assay
Luciferase Solution (Green cap)	10 µl×1	
ATP Assay Buffer	5.5 ml×1	
Oligomycin (Yellow cap)	×1	

Preparation of Oligomycin solution

(1) Dilute the 10 mmol/l Oligomycin stock solution 1,000-fold with culture medium (to give 10 μmol/l Oligomycin solution).

※ For example, to prepare 1 ml of Oligomycin solution, mix 1 μl of 10 mmol/l Oligomycin stock solution and 1 ml of culture medium.

(2) Dilute the solution prepared in step (1) 4-fold with culture medium to give 2.5 μmol/l Oligomycin solution.

(3) Dilute the solution prepared in step (2) 2-fold with culture medium to give 1.25 μmol/l Oligomycin solution.

※ Where cells are to be stimulated with a drug, add the drug to the medium used to prepare the Oligomycin solution so that the drug concentration is the same as required for stimulation (see Figure 5).

※ Perform a preliminary experiment and determine whether to use 2.5 or 1.25 μmol/l Oligomycin solution in the experiment (see Experimental Procedure-Preliminary Experiment (optimization of cell number and Oligomycin concentration)).

※ Use Oligomycin solution within the same day (it cannot be stored).

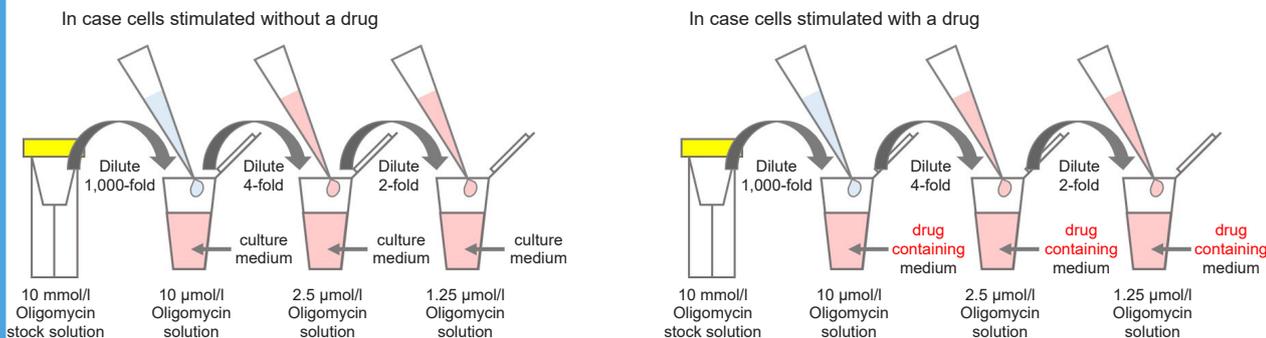
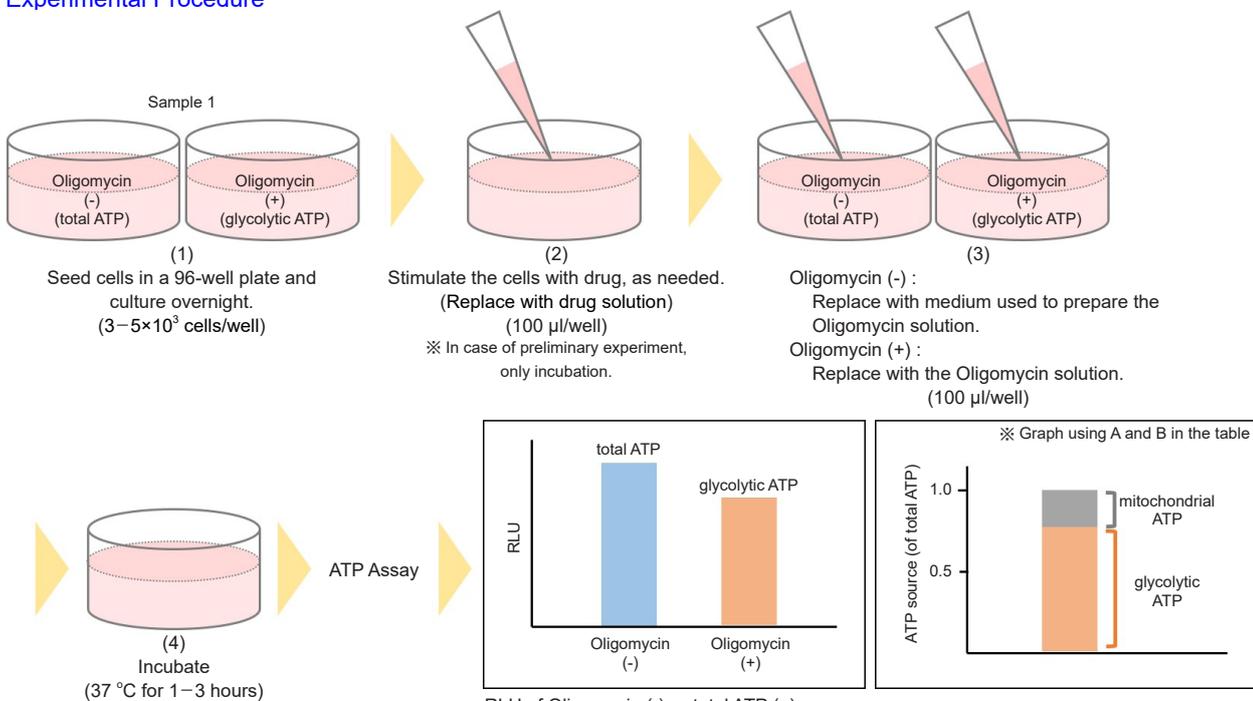


Figure 5. Preparation of Oligomycin-containing medium

Experimental Procedure



RLU of Oligomycin (-) = total ATP (a)
 RLU of Oligomycin (+) = glycolytic ATP (b)
 (RLU of Oligomycin (-)) - (RLU of Oligomycin (+)) = mitochondrial ATP (c)

total ATP (a)	(a)/(a) = 1	
glycolytic ATP (b)	(b)/(a)	A
mitochondrial ATP(c)	1 - {(b)/(a)}	B

— Preliminary Experiment (optimization of cell number and Oligomycin concentration) —

(1) Seed cells in a 96-well white microplate (3–5×10³ cells/well) and culture in an incubator (37 °C, 5% CO₂).

※ Adjust the optimal cell number to be seeded according to the drug stimulation time (see Table 5).

※ See Figure 6 for an assay plate format when performing the optimization.

※ Use the same incubation time as the time required for drug stimulation.

For example, if the drug stimulation time is 24 hours, culture cells overnight after seeding, and then culture for another 24 hours.

(2) Remove the medium and add 100 μl of Oligomycin solution to test sample Oligomycin (+) wells.

(3) Add 100 μl of the medium used to prepare the Oligomycin solution to test sample Oligomycin (-) wells.

(4) Incubate the microplate for 1–3 hours in an incubator (37 °C, 5% CO₂).

※ For incubation time 1–3 hours, the results are almost the same.

(5) Perform the ATP Assay.

(6) Compare the luminescence value of with and without Oligomycin treatment.

※ Choose the optimum cell number and Oligomycin concentration by referring to Table 6 and Figure 7.

Table 5. Example of cell number

	Drug treatment time	
	without	24 hours
HeLa, HepG2, A549	5–10×10 ³ cells	3–5×10 ³ cells
WI-38	5–10×10 ¹⁰ cells	5×10 ³ cells

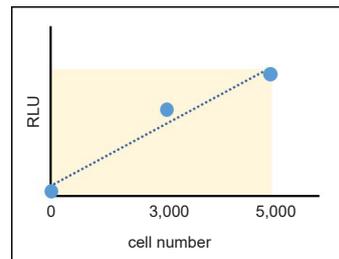
	1	2	3
A	Oligomycin (-) 5,000 cells/well		
B	Oligomycin (-) 3,000 cells/well		
C	blank (medium only)		
D	2.5 $\mu\text{mol/l}$ Oligomycin 5,000 cells/well		
E	1.25 $\mu\text{mol/l}$ Oligomycin 5,000 cells/well		
F	2.5 $\mu\text{mol/l}$ Oligomycin 3,000 cells/well		
G	1.25 $\mu\text{mol/l}$ Oligomycin 3,000 cells/well		

: For optimization of cell number
 : For optimization of Oligomycin concentration (5,000 cells/well)
 : For optimization of Oligomycin concentration (3,000 cells/well)

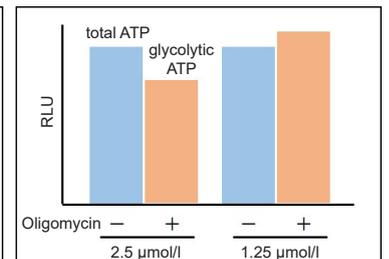
Figure 6.
Example of an assay plate format for optimization of cell number and Oligomycin concentration (n=3)

Table 6. Example of Oligomycin concentration

	Oligomycin treatment time	
	1 hour	3 hours
HepG2, A549, WI-38	1.25 or 2.5 $\mu\text{mol/l}$	1.25 or 2.5 $\mu\text{mol/l}$
HeLa	2.5 $\mu\text{mol/l}$	1.25 or 2.5 $\mu\text{mol/l}$



Choose a suitable cell number from within the linear portion of the plotted curve.



Choose a suitable Oligomycin concentration which the RLU of glycolytic ATP does not exceed the RLU of total ATP as much as possible.

Figure 7.
Determination of suitable cell number and Oligomycin concentration

— Experiment —

- Seed cells in a 96-well white microplate (choose a suitable cell number by optimization) and culture overnight in an incubator (37 °C, 5% CO₂).
※ See Figure 8 for an assay plate format.
- Remove the medium and stimulate the cells with the drug, as needed.
※ Add drug solution to wells.
- Remove the medium and add 100 μl of Oligomycin solution (choose a suitable Oligomycin concentration by optimization) to test sample Oligomycin (+) wells.
- Add 100 μl of the medium used to prepare the Oligomycin solution to test sample Oligomycin (-) wells.
- Incubate the microplate for 1–3 hours in an incubator (37 °C, 5% CO₂).
※ For incubation time 1–3 hours, the results are almost the same.
- Perform the ATP Assay.
- Compare the luminescence value of with and without Oligomycin treatment.

	1	2	3
A	Sample 1 Oligomycin (-)		
B	Sample 1 Oligomycin (+)		
C	Sample 2 Oligomycin (-)		
D	Sample 2 Oligomycin (+)		

Figure 8.
Example of an assay plate format (n=3)

Experimental Example : Metabolism shift of glycolytic inhibitor-treated HeLa cells

- HeLa cells were seeded in a 96-well white microplate (5×10³ cells/well in MEM containing 10% fetal bovine serum and 1% penicillin-streptomycin) and cultured overnight in an incubator (37 °C, 5% CO₂).
※ Six wells were prepared for glycolytic inhibitor stimulation and six wells for the control treatment.
- The medium was removed, and glycolytic inhibitor solution (100 μl , in MEM containing 10% FBS) was added to six cell-seeded wells of the 96-well microplate, and MEM containing 10% FBS (100 μl) was added to the other six cell-seeded wells.
- The cells were cultured overnight in an incubator (37 °C, 5% CO₂).
- The medium was removed, and 1.25 $\mu\text{mol/l}$ Oligomycin solution (100 μl) was added to the Oligomycin (+) wells.
- The medium (100 μl) used to prepare the Oligomycin solution was added to the Oligomycin (-) wells.
- The cells were cultured for 3 hours in an incubator (37 °C, 5% CO₂).
- After incubation, ATP working solution (100 μl) was added to each well.
- The assay plate was orbitally shaken for 2 minutes in a microplate reader.
- The assay plate was incubated at 25 °C for 10 minutes in a microplate reader.
- Luminescence was measured using a plate reader, and total ATP, glycolytic ATP and mitochondrial ATP were calculated.

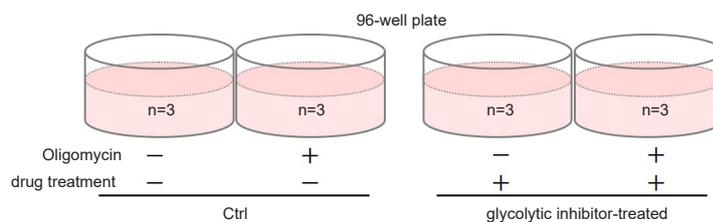
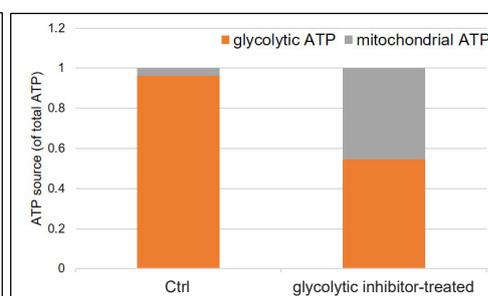
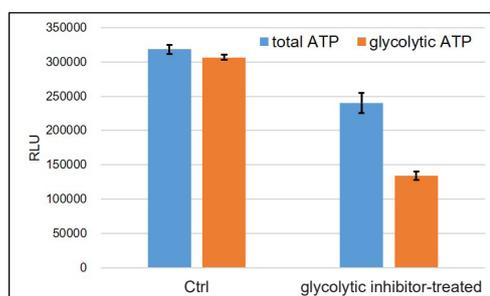
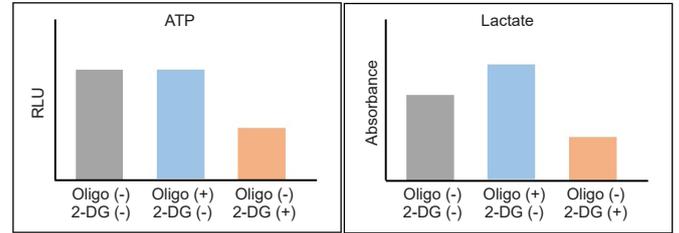
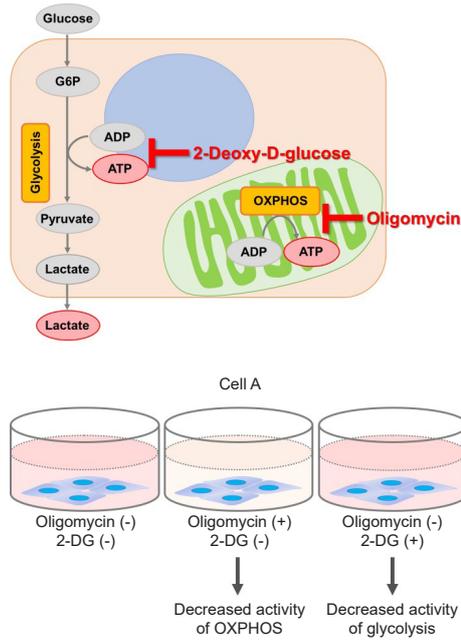


Figure 9.
Metabolism shift of glycolytic inhibitor-treated HeLa cells
Glycolytic inhibitor treatment caused a metabolism shift in HeLa cells from glycolysis to oxidative phosphorylation.



Overview of Analysis

This is a protocol to know whether the cells are mainly dependent on glycolysis or oxidative phosphorylation by measuring ATP and lactate production. Oligomycin and 2-deoxy-D-glucose (2-DG) inhibits ATP synthesis in oxidative phosphorylation and glycolysis, respectively.



ATP : Decrease by glycolytic inhibition
: No change by OXPHOS inhibition

Lactate : Decrease by glycolytic inhibition
: Increase by OXPHOS inhibition

From the change in lactate, the glycolysis is activated when oxidative phosphorylation is inhibited. Furthermore, although ATP decrease when glycolysis is inhibited, ATP does not change even when oxidative phosphorylation is inhibited. So, these are suggested that the cells are mainly dependent on glycolysis.

Reagents

Table 7. Reagents for analysis of metabolic pathway dependence

Dye Mixture (Blue cap)	×1	Use for Lactate Assay
Lactate Standard (10 mmol/l) (Blue cap)	150 μl×1	
LDH Solution (Purple cap)	12 μl×1	
Lactate Assay Buffer	5.5 ml×1	
Reconstitution Buffer (White cap)	550 μl×1	
Substrate (Green cap)	×1	Use for ATP Assay
Luciferase Solution (Green cap)	10 μl×1	
ATP Assay Buffer	5.5 ml×1	
Oligomycin (Yellow cap)	×1	Use as an inhibitor
2-DG (Red cap)	140 μl×1	

Preparation of Oligomycin solution

(1) Dilute the 10 mmol/l Oligomycin stock solution 1,000-fold with culture medium (to give 10 μmol/l Oligomycin solution).

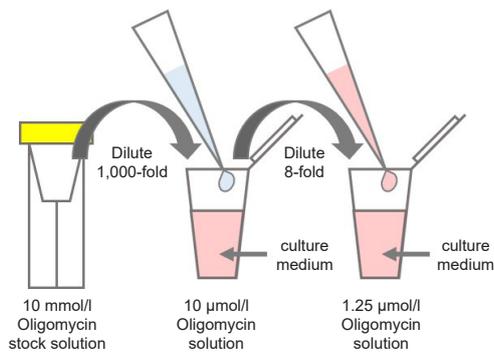
※ For example, to prepare 1 ml of Oligomycin solution, mix 1 μl of 10 mmol/l Oligomycin stock solution and 1 ml of culture medium.

(2) Dilute the solution prepared in step (1) 8-fold with culture medium to give 1.25 μmol/l Oligomycin solution.

※ Where cells are to be stimulated with a drug, add the drug to the medium used to prepare the Oligomycin solution so that the drug concentration is the same as required for stimulation (see Figure 10).

※ Use Oligomycin solution within the same day (it cannot be stored).

In case cells stimulated without a drug



In case cells stimulated with a drug

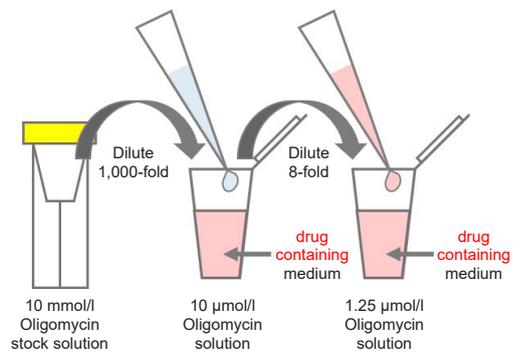


Figure 10. Preparation of Oligomycin-containing medium

Preparation of 2-DG solution

Dilute the 2-DG (red cap) 40-fold with culture medium (to give 22.5 mmol/l 2-DG solution).

- ※ For example, to prepare 1 ml of 2-DG solution, mix 25 μ l of 2-DG and 975 μ l of culture medium.
- ※ Where cells are to be stimulated with a drug, add the drug to the medium used to prepare the 2-DG solution so that the drug concentration is the same as required for stimulation (see Figure 11).
- ※ Use 2-DG solution within the same day (it cannot be stored).

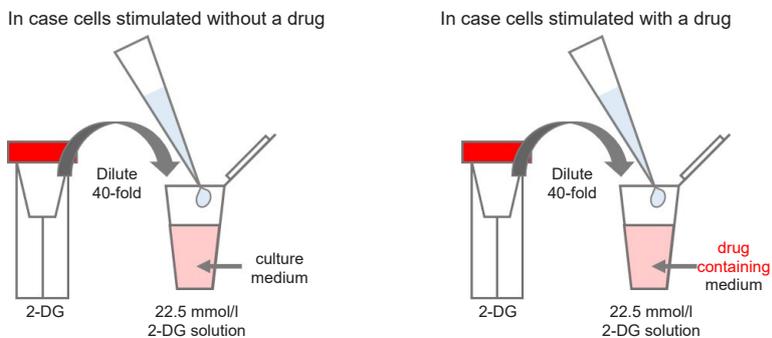
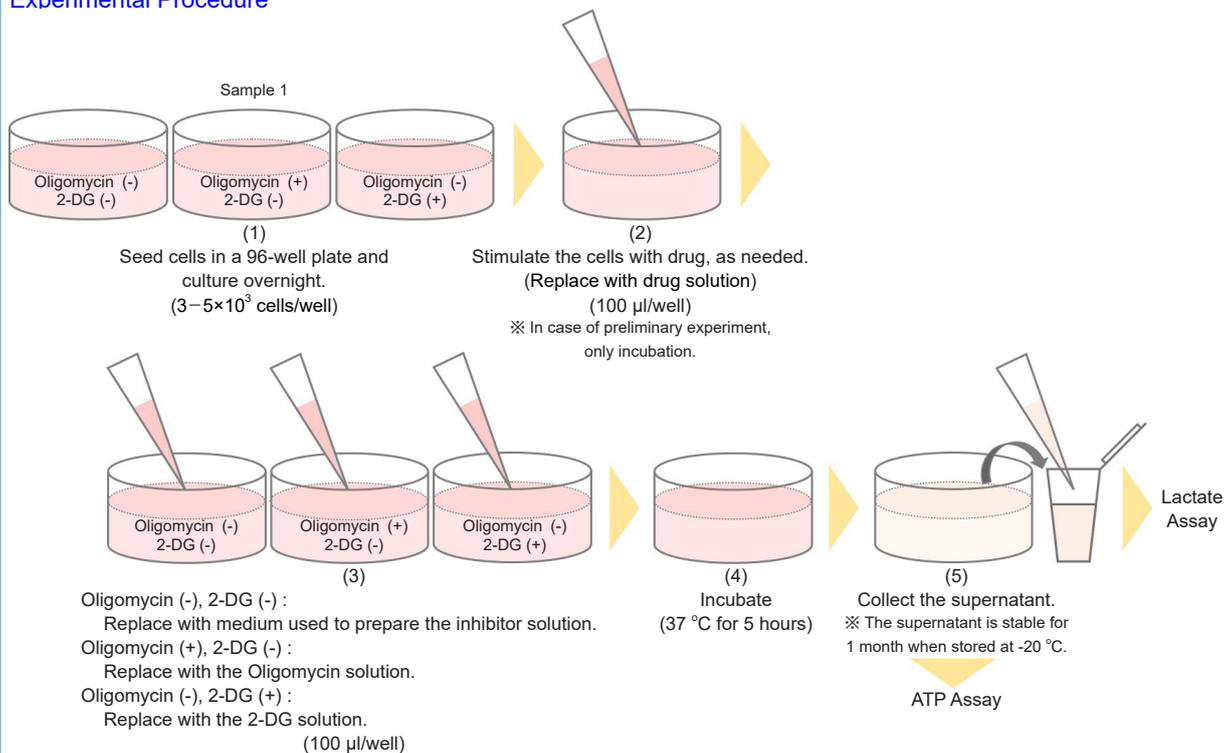


Figure 11. Preparation of 2-DG-containing medium

Experimental Procedure



— Preliminary Experiment (optimization of cell number) —

- Seed cells in a 96-well white microplate (3–5×10³ cells/well) and culture in an incubator (37 °C, 5% CO₂).
 ※ Adjust the optimal cell number to be seeded according to the drug stimulation time (see Table 8).
 ※ See Figure 12 for an assay plate format when performing the optimization.
- Use the same incubation time as the time required for drug stimulation.
 For example, if the drug stimulation time is 24 hours, culture cells overnight after seeding, and then culture for another 24 hours.
- Remove the medium and add 100 μ l of fresh medium.
- Incubate the microplate for 5 hours in an incubator (37 °C, 5% CO₂).
- Perform the ATP Assay.
- Compare the luminescence value and choose the suitable cell number (see Figure 13).

	1	2	3
A	5,000 cells/well		
B	3,000 cells/well		
C	blank (medium only)		

Figure 12.

Example of an assay plate format for optimization of cell number (n=3)

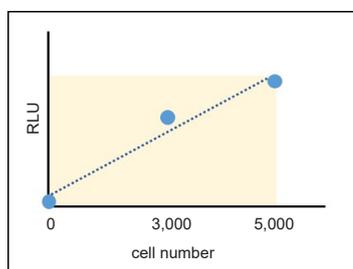


Table 8. Example of cell number

	Drug treatment time	
	without	24 hours
HeLa, HepG2, A549	5–10×10 ³ cells	3–5×10 ³ cells
WI-38	5–10×10 ³ cells	5×10 ³ cells

Figure 13.

Determination of suitable cell number
 Choose a suitable cell number from within the linear portion of the plotted curve.

— Experiment —

- (1) Seed cells in a 96-well white microplate (choose a suitable cell number by optimization) and culture in an incubator (37 °C, 5% CO₂).
- ※ See Figure 14 for an assay plate format.
- (2) Remove the medium and stimulate the cells with the drug, as needed.
- ※ Add drug solution to wells.
- (3) Remove the medium and add 100 µl of 1.25 µmol/l Oligomycin solution to test sample Oligomycin (+), 2-DG (-) wells.
- (4) Add 100 µl of 2-DG solution to test sample Oligomycin (-), 2-DG (+) wells.
- (5) Add 100 µl of the medium used to prepare the Oligomycin solution and 2-DG solutions to test sample Oligomycin (-), 2-DG (-) wells.
- (6) Incubate the microplate for 5 hours in an incubator (37 °C, 5% CO₂).
- (7) Collect 20 µl of each cell culture supernatant into a 1.5 ml microtube and perform the Lactate Assay.
- ※ See Figure 15 for an assay plate format when performing the Lactate Assay.
- (8) Perform the ATP Assay using the cells remaining in the 96-well white microplate.
- (9) Compare the luminescence and absorbance values of with and without Oligomycin and 2-DG treatment, respectively.

	1	2	3
A	Sample 1 Oligomycin (-), 2-DG (-)		
B	Sample 1 Oligomycin (+), 2-DG (-)		
C	Sample 1 Oligomycin (-), 2-DG (+)		
D	Sample 2 Oligomycin (-), 2-DG (-)		
E	Sample 2 Oligomycin (+), 2-DG (-)		
F	Sample 2 Oligomycin (-), 2-DG (+)		

Figure 14.

Example of an assay plate format for seeding cells
(n=3)

	1	2	3
A	Sample 1 Oligomycin (-), 2-DG (-)		
B	Sample 1 Oligomycin (+), 2-DG (-)		
C	Sample 1 Oligomycin (-), 2-DG (+)		
D	Sample 2 Oligomycin (-), 2-DG (-)		
E	Sample 2 Oligomycin (+), 2-DG (-)		
F	Sample 2 Oligomycin (-), 2-DG (+)		
G	medium only (background)		

Figure 15.

Example of an assay plate format for Lactate Assay
(n=3)

Experimental Example : Metabolic pathway dependence of glycolytic inhibitor-treated HeLa cells

- (1) HeLa cells were seeded in a 96-well white microplate (3×10³ cells/well in MEM containing 10% fetal bovine serum and 1% penicillin-streptomycin) and cultured overnight in an incubator (37 °C, 5% CO₂).
- ※ Nine wells were prepared for glycolytic inhibitor stimulation and nine wells for the control treatment.
- (2) The medium was removed, and glycolytic inhibitor solution (100 µl, in MEM containing 10% FBS) was added to nine wells of the 96-well microplate, and MEM containing 10% FBS (100 µl) was added to the other nine wells.
- (3) The cells were cultured overnight in an incubator (37 °C, 5% CO₂).
- (4) The medium was removed, and 1.25 µmol/l Oligomycin solution (100 µl) was added to Oligomycin (+), 2-DG (-) wells.
- (5) The 22.5 mmol/l 2-DG solution (100 µl) was added to Oligomycin (-), 2-DG (+) wells.
- (6) The medium (100 µl) used to prepare the Oligomycin and 2-DG solutions was added to Oligomycin (-), 2-DG (-) wells.
- (7) The cells were cultured for 5 hours in an incubator (37 °C, 5% CO₂).
- (8) After incubation, 20 µl of each cell culture supernatant was transferred to a 1.5 ml microtube and diluted 10-fold with ddH₂O to prepare sample solutions for the Lactate Assay.
- ※ For background control, MEM only was also diluted 10-fold with ddH₂O.
- (9) ATP working solution (100 µl) was added to the cells remaining in each well of the 96-well white microplate.
- (10) The assay plate was orbitally shaken for 2 minutes in a microplate reader.
- (11) The assay plate was incubated at 25 °C for 10 minutes in a microplate reader.
- (12) Luminescence was measured using a plate reader, and the luminescence values of with and without Oligomycin and 2-DG treatments were compared.
- (13) The sample solutions (20 µl each) prepared in step (8) were added to individual wells of a 96-well microplate.
- (14) Lactate working solution (80 µl) was added to each well.
- (15) The assay plate was incubated at 37 °C for 30 minutes.
- (16) The absorbance at 450 nm was measured using a microplate reader, and the absorbance values of with and without Oligomycin and 2-DG treatments were compared.

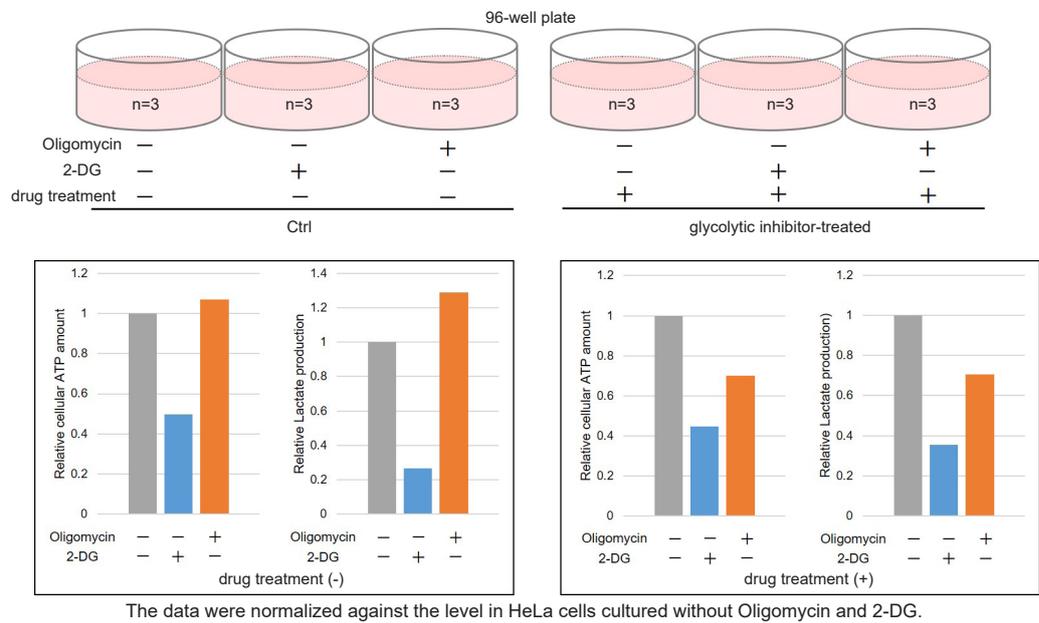


Figure 16.

Comparison of metabolic pathway dependence of glycolytic inhibitor-treated and non-treated (Ctrl) HeLa cells

Without glycolytic inhibitor-treatment, the metabolism of HeLa cells is mainly dependent on glycolysis, but glycolytic inhibitor-treated HeLa cells are dependent on oxidative phosphorylation.

References

- 1) R. Shiratori, et al., *Sci. Rep.*, **2019**, *9*, 18699.
- 2) J. Kim, et al., *Cancer Cell*, **2019**, *35*, 191-203.
- 3) H. Kobayashi, et al., *Nature. Chem. Bio.*, **2021**, *17*, 335-343.

For research use only. Not for use in diagnostic procedures.
If you need more information, please contact Dojindo technical service.

Dojindo Laboratories
2025-5 Tabaru, Mashiki-machi, Kamimashiki-gun, Kumamoto
861-2202, Japan Phone: +81-96-286-1515 Fax: +81-96-286-1525
E-mail: info@dojindo.co.jp Web: www.dojindo.co.jp

Dojindo Molecular Technologies, Inc.
Tel: +1-301-987-2667 Web: <http://www.dojindo.com/>
Dojindo EU GmbH
Tel: +49-89-3540-4805 Web: <http://www.dojindo.eu.com/>
Dojindo China Co., Ltd
Tel: +86-21-6427-2302 Web: <http://www.dojindo.cn/>
Tel: +86-10-8225-1765

G270 : Glycolysis/OXPHOS Assay Kit