



FCCP Optimization using the Cell Mito Stress Test with the Agilent Seahorse XF^e96/XF96

Basic Procedure

NOTE: For this assay, seed cells at the optimal cell number.

The Cell Mito Stress Test is run with six different concentrations of FCCP to determine the optimal FCCP concentration to use in your XF assays.

Plate Layout

[FCCP] 0 μ M 0.125 μ M 0.25 μ M 0.5 μ M 1.0 μ M 2.0 μ M

	1	2	3	4	5	6	7	8	9	10	11	12
A	Black	Green	Purple	Purple	Yellow	Yellow	Red	Red	Cyan	Cyan	Magenta	Black
B	Green	Green	Purple	Purple	Yellow	Yellow	Red	Red	Cyan	Cyan	Magenta	Magenta
C	Green	Green	Purple	Purple	Yellow	Yellow	Red	Red	Cyan	Cyan	Magenta	Magenta
D	Green	Green	Purple	Purple	Yellow	Yellow	Red	Red	Cyan	Cyan	Magenta	Magenta
E	Green	Green	Purple	Purple	Yellow	Yellow	Red	Red	Cyan	Cyan	Magenta	Magenta
F	Green	Green	Purple	Purple	Yellow	Yellow	Red	Red	Cyan	Cyan	Magenta	Magenta
G	Green	Green	Purple	Purple	Yellow	Yellow	Red	Red	Cyan	Cyan	Magenta	Magenta
H	Black	Green	Purple	Purple	Yellow	Yellow	Red	Red	Cyan	Cyan	Magenta	Black

Injections

NOTE: All compound injection solutions are to be made at 10x their final respective concentrations in the wells.



Agilent Technologies

Port A: oligomycin

Port A: oligomycin: 1.0 μM (optimal) final concentration in the well

Port B: FCCP

1. Columns 1-2: 0 μM final concentration in the well
2. Columns 3-4: 0.125 μM final concentration in the well
3. Columns 5-6: 0.25 μM final concentration in the well
4. Columns 7-8: 0.50 μM final concentration in the well
5. Columns 9-10: 1.0 μM final concentration in the well
6. Columns 11-12: 2.0 μM final concentration in the well

Port C: rotenone/antimycin A: 0.5 μM final concentration in the well

Protocol

1. Warm the Cell Mito Stress Test Assay Medium to 37°C. For information on assay media preparation, see: http://www.agilent.com/cs/library/usermanuals/public/XFe96_DAY_OF_MEDIA_PREP.pdf and http://www.agilent.com/cs/library/usermanuals/public/XF96_DAY_OF_MEDIA_PREP.pdf. Adjust pH to 7.4 ± 0.1 at 37°C.
2. Retrieve the cell plate from the CO₂ incubator.
3. Look at cells under the microscope to:
 - a. Confirm cell health, morphology, seeding uniformity and purity (no contamination)
 - b. Ensure cells are adhered, and show a consistent monolayer.
 - c. Make sure no cells were plated in the background correction wells.
4. Wash cells two times with Cell Mito Stress Test Assay Medium, for detailed instructions see: http://www.agilent.com/cs/library/usermanuals/public/XFe96_DAY_OF_WASHING_CELLS.pdf or http://www.agilent.com/cs/library/usermanuals/public/XF96_DAY_OF_WASHING_CELLS.pdf.
5. Look at cells under the microscope to ensure that cells were not disturbed or washed away.
6. Place the plate in a 37°C incubator without CO₂ for 45-60 min prior to the assay.
7. Prepare Stock Compounds for Injection.
 - a. Important: Use compounds the same day they are reconstituted. Do not refreeze. Discard any remaining compound.
 - b. Remove one foil pouch from XF Cell Mito Stress Test Kit box.
 - c. Remove the three tubes containing oligomycin (blue cap), FCCP (yellow cap), and rotenone/antimycin A (red cap).



Agilent Technologies

- d. Resuspend contents of each tube with assay medium using volumes described in the table below with a p1000 pipette to make Stock solutions for each compound. Gently pipette up and down (~10 times) or vortex with cap on to solubilize the compounds.

	Volume of Assay Medium	Final Stock Concentration
Oligomycin	630 μ L	100 μ M
FCCP	720 μ L	100 μ M
Rotenone/AntimycinA	540 μ L	50 μ M

8. Prepare the compound solutions that will be loaded into the cartridge injection ports.
a. Prepare 3 mL of 10 μ M oligomycin in assay medium:

Port A Oligomycin	[Final well] (μ M)	Stock volume (μ L)	Medium volume (μ L)
	1.0	300	2,700

- b. Prepare serial dilutions of FCCP in assay medium, starting with Tube a as detailed below:

Port B FCCP	Tube	[Final well] (μ M)	Stock volume (μ L)	Medium volume (μ L)
	a	2.0	700	2,700
	b	1.0	1500 from a	1500
	c	0.5	1500 from b	1500
	d	0.25	1500 from c	1500
	e	0.125	1500 from d	1000
	f	0	0	2000

- c. Pipette 300 μ L of the rotenone/antimycin A stock into a 2700 μ L aliquot of assay medium:

Port C Rotenone/Antimycin A	[Final well] (μ M)	Stock volume (μ L)	Medium volume (μ L)
	0.5	300	2,700

9. Obtain a hydrated cartridge from the non-CO₂ incubator. Load the cartridge in each port as outlined below.
a. Port A – 1.0 μ M oligomycin final concentration in the well. **Load 20 μ L** of the 10x stock into each Port A.
b. Port B – FCCP dilutions: Note the layout! **Load 22 μ L** of each 10x solution into the B ports in the appropriate columns shown below.
i. Columns 1-2: 0 μ M final concentration in the well
ii. Columns 3-4: 0.125 μ M final concentration in the well
iii. Columns 5-6: 0.25 μ M final concentration in the well



- iv. Columns 7-8: 0.50 μM final concentration in the well
 - v. Columns 9-10: 1.0 μM final concentration in the well
 - vi. Columns 11-12: 2.0 μM final concentration in the well
- c. Port C – 0.5 μM Rot/AA final concentration in the well. **Load 20 μL** of the 10x stock into each Port C.
10. Create or load your assay design on the XF[®]96/XF96 Controller. Default Mix-Wait-Measure times are 3 min – 2 min – 3 min. In general, 3 basal rate measurements are taken prior to the first injection; then 3 rate measurements after each injection.
 11. When ready, initiate the cartridge calibration and follow the screen prompts and load the cartridge.
 12. After cartridge calibration is complete, follow the software prompts to replace the Utility Plate with the cell culture and initiate the assay.

Learn more

www.agilent.com/en-us/promotions/seahorse-xf-technology

For Research Use Only. Not for use in diagnostic procedures.

Product specifications and descriptions in this document are subject to change without notice.

© Agilent Technologies, Inc., 2014, 2017

Published in the USA, June 30, 2017



Agilent Technologies