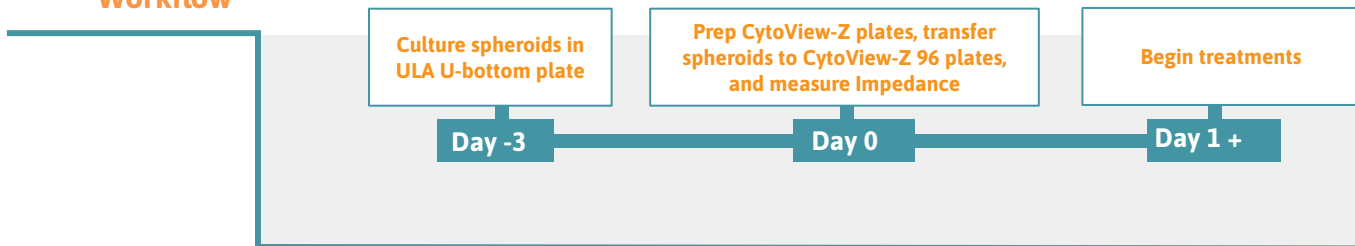


Cell Culture Protocol

Cancer Spheroids for the Maestro Z

Workflow



Culturing Spheroids

1. Thaw and culture cancer cells in accordance with supplier recommendations and passage as needed.
2. Remove flask(s) of cells from the incubator, aspirate the media, and rinse with PBS. Using trypsin, or another cell dissociating agent, detach and collect the cells from the flasks per reagent recommendations.
3. Remove a sample of the cell suspension and count the cells using a hemocytometer to determine the total number of viable cells.
4. Transfer the cell suspension to a 15 mL conical tube and centrifuge to a pellet. Aspirate the supernatant, being careful not to disturb the cell pellet.
5. Dilute the cell suspension in complete medium to a working concentration of cells per 200 μL .
6. Transfer 200 μL of the cell suspension to each well of a ultra-low attachment (ULA) U-bottom plate to achieve the desired number of cells per spheroid.
7. Centrifuge the plate at 200 x g for 5 minutes then place in the incubator at 37°C and 5% CO_2 .
8. Allow the spheroids to grow for 4 days.

Tip

- For non-ULA plates:
1. Add 200 μL of STEMCell Technologies™ Anti-Adherence Rinsing Solution to each well.
 2. Centrifuge at 200 x g for 5 minutes.
 3. Aspirate and wash with PBS.
 4. Add 200 μL of PBS to each well.
 5. Aspirate the PBS when the cell suspension is ready to transfer.

Tip

Use "mid-low" speed for centrifuge deceleration to avoid cell spreading against the walls of the wells.

Preparation of CytoView-Z Plate and Baseline Recording

9. After 4 days of spheroid culture, add 100 μL of extracellular matrix (ECM) to the CytoView-Z 96 plate.
10. Incubate the plate at 37°C and 5% CO_2 for at least 1 hour.
11. Aspirate the excess ECM from the plate.
12. Add 100 μL of complete medium to the wells of the plate and 8 mL of sterile water to the on-plate reservoirs to increase humidity.
13. Dock the plate in a Maestro platform to measure the media-only baseline. Transfer the plate to a biosafety cabinet when the baseline is complete.

Tip

Reconstitute ECM as recommended by supplier. Store as single use aliquots for more consistent results.



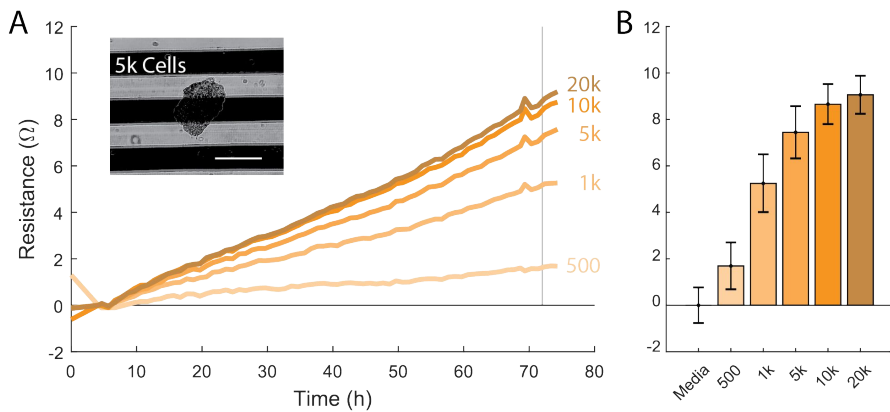


Figure 1: Monitoring attachment and proliferation of cancer spheroids

(A) A549 spheroids plated on a CytoView-Z 96-well plate from low to high density (500 to 20k cells/spheroid). Inset: A549 spheroid formed from 5k cells. Scale bar = 200 μ m. (B) Resistance at 72 hours post-plating for A549 spheroids.

Transfer Spheroids to CytoView-Z Plate

14. Check the spheroids under a microscope for stability and cell-to-cell adhesion. The spheroids should have a defined shape (e.g., no single cell scatter).
15. Insert a P-1000 micropipette with wide bore tips into the U-bottom well and make sure the tip is touching the bottom of the well. The tip should surround the spheroid.
16. Pipette up and down a few times to loosen the spheroid from the U-bottom well.
17. With the tip still touching the bottom, pipette up 100 μ L from the U-bottom well and transfer the media with the spheroid into the well of the CytoView-Z plate.

Note: For multiple spheroids per well, combine spheroids into a single U-bottom well. Wait for each spheroid to settle to the bottom. Then, pipette all the spheroids in 100 μ L and transfer to the CytoView-Z plate.

18. Dock the plate into the Maestro platform.

Tip

Periodically check if the transfer was successful. If there is no spheroid in the CytoView-Z well:

1. Check the coordinating U-bottom well for the spheroid(s).
2. Repeat steps 15-17.
3. Re-check the U-bottom and CytoView-Z wells under a microscope.

Tip

CytoView-Z wells include a viewing window. If spheroid(s) land in the area without electrodes (dark bands), it is advised to disable that well in Axis Z.

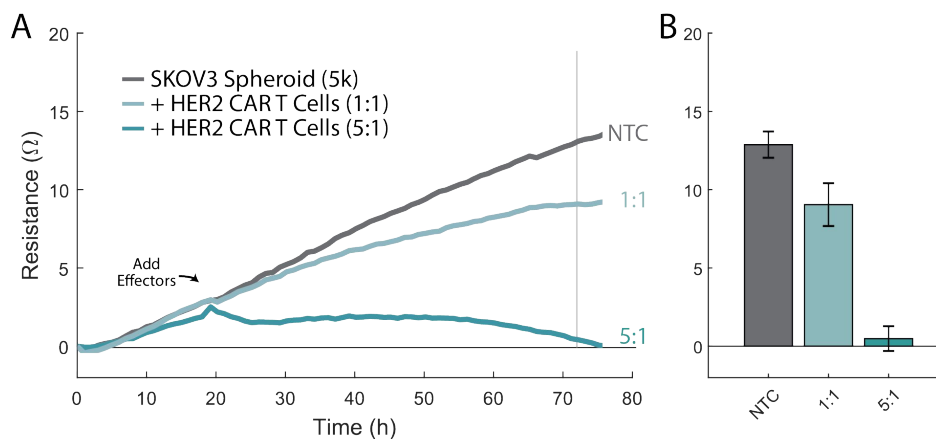


Figure 2: Immune cell-mediated killing of SKOV3 spheroids

(A) SKOV3 spheroids (5k cells per spheroid) were plated on a CytoView-Z 96 plate. HER2-specific CAR T cells were added at 1:1 and 5:1 effector-to-target cell ratios after 20 hours. (B) Resistance at 52 hours post-effector addition.

Required Materials

Consumables

Item	Vendor
CytoView-Z 96 Plate	Axion BioSystems
Extracellular matrix	Various
Trypsin-EDTA (0.5%) or other dissociation reagent	Various
Gibco™ PBS, pH 7.4 Ca and Mg-free	Various
1000 μ L Wide Bore Pipette Tips	Thomas Scientific (#1141M27)
ULA U-Bottom Plates	Corning (#4515)
Cell Culture Media	Various
15 mL Conical Tubes	Various

Equipment

Item	Vendor
Maestro Edge, Pro, Z, or ZHT	Axion BioSystems
Biological Safety Cabinet	Various
Cell Culture Incubator	Various
P1000 micropipette	Various
Hemocytometer or Cell Counter	Various
Phase Contrast Microscope	Various
Tabletop Centrifuge	Various
37°C Water Bath	Various