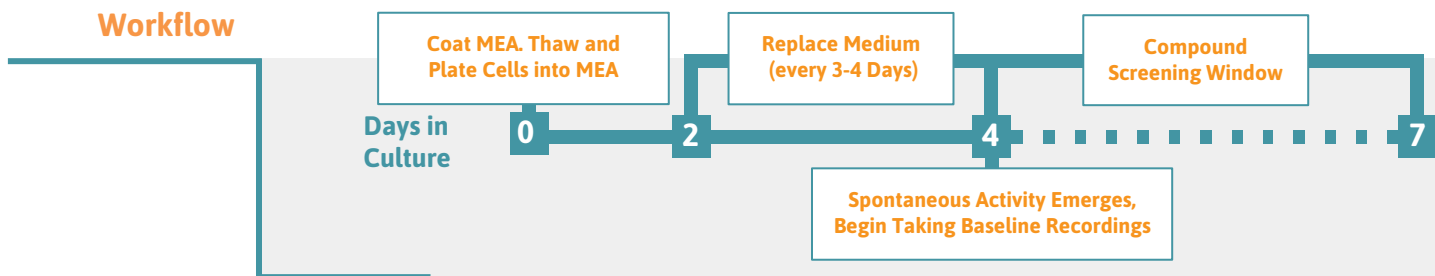


Cell Culture Protocol

QBM Cell Science Rat Dorsal Root Ganglion Neurons



Preparing the MEA Plate

1. Place a 5 μ l droplet of 0.1 % PEI solution over the recording electrode area of each well in the MEA plate. See Figure 1 on page 2 for appropriate drop placement.
2. Incubate the PEI-coated MEA plate in a cell culture incubator at 37°C, 5% CO₂ for at least 60 minutes.
3. Rinse PEI from the culture surface with 200 μ L of sterile DI water 4 times, then allow the MEA plate to air dry overnight.

Tip

Recommended to add 6-8 mL of sterile water to the on-plate reservoirs to increase humidity.

Culturing QBM Rat Cortical Neurons

4. Thaw Rat DRG Neurons by holding the tube partially submerged in a 37°C water bath for 3 minutes, making sure to not submerge the cap. Spray the outside with 70% ethanol and transfer to a biosafety cabinet. Dilute the cells with 37°C complete media.
5. Remove a sample of the cell suspension and count the neurons using a hemocytometer to determine both the viability and total number of viable cells. Transfer the cell suspension to a 15 ml conical tube.
6. Centrifuge the cell suspension at 100 x g for 1 minute.
7. Aspirate the supernatant, being careful not to disturb the cell pellet.
8. Dilute the cell suspension in Complete PNGM™ BulletKit™ Neuron Growth Medium + NGF- β and laminin (20 μ g/mL). The final cell concentration should be 16,000,000 viable neurons/ml.

Tip

Ensure the neurons are evenly suspended before removing an aliquot to count.

Tip

The rat DRG neurons are sensitive to centrifugation, so care should be taken to monitor speed and duration during this step. The cell provider does not recommend centrifugation and is not responsible for cell death induced by centrifugation.

Plating QBM Rat Cortical Neurons onto the MEA

9. Place a 5 μ l droplet of DRG Neuron suspension over the recording electrode area of each well of the MEA. See Figure 1 on page 2 for appropriate drop placement.
10. Incubate the MEA plate with the seeded neurons in a cell culture incubator at 37°C, 5% CO₂ for 1 hour.

Tip

Prepare the laminin fresh from frozen aliquots for every cell culture.

11. Gently add 1/2 of the final volume of culture medium to each well of the MEA. Adding the medium too quickly will dislodge the adhered neurons. Recommended well volumes for each plate type are: 6- and 12-well = 1000 μ l, 24-well = 500 μ l, 48-well = 300 μ l, 96 well = 200 μ l.
12. Repeat step 11 a second time to reach the final recommended volume of medium.
13. Incubate in a cell culture incubator at 37°C, 5% CO₂.
14. *Optional:* To control Schwann cell population, add antimetabolites 4 hours after seeding into medium, 17.5 μ g/mL uridine and 7.5 μ g/mL of 5-fluoro-2'-deoxyuridine. Continue the antimetabolite addition to the complete medium with each maintenance medium change.
15. For optimal cell health, be sure to exchange 50% of the medium every 3-4 days. Though neural spikes may be detectable within 2 days, optimal neural activity is typically achieved after 4 days in culture.

Tip

Using a pipette, add medium first in a semi-circle along the outer edge of the well. Progressively add medium to either side of the well so it fills evenly towards the center. The goal is to prevent a rush of medium in either direction that might dislodge the neurons.

Drop Placement

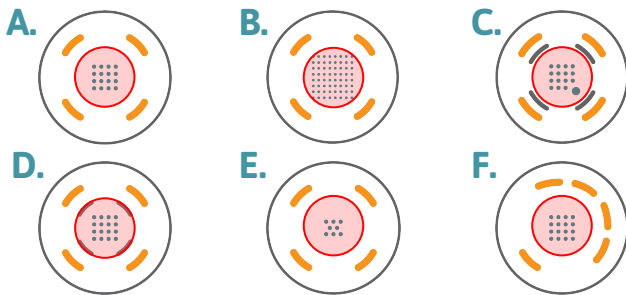


Figure 1: Drop Placement Diagram

The layouts above represent the bottom surfaces of wells in (A) a 48-well MEA, (B) a 6- or 12-well MEA, (C) a 24-well MEA or 48 well E-Stim+ MEA, (D) a 48-well AccuSpot MEA, (E) a 96-well MEA, and (F) a 48-well CytoView MEA. The number of electrodes per well is different across the plate formats, however the drop placement is the same, with the drop (red circle) centered on the recording electrodes and staying within the ground electrodes. On plate types (C) with the addition of the stim-paddle in the lower right corner of the array, it is important to make sure the droplet covers this feature. The droplet may need to be manipulated after placement of the pre-treatment to ensure stim-paddle coverage.

Visualization of Typical Neuron Seeding Results

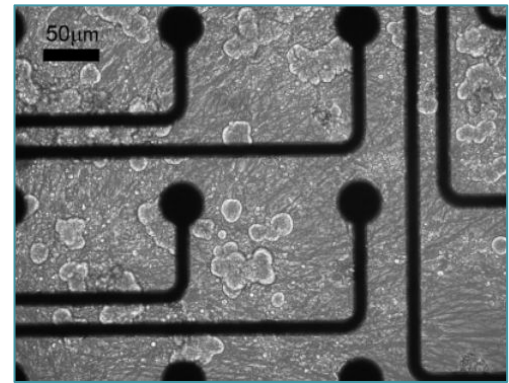


Figure 2: QBM Rat DRG Neuron Morphology

Primary Rat DRG Neurons (80,000) at day 6 *in vitro* in a 12-well CytoView MEA, 20x magnification. Notice that neuron morphology is easily recognizable.

Required Materials

Consumables

Item	Vendor	Catalog #
Axion MEA (6, 12, 24, 48, or 96-Well)	Axion BioSystems	
QBM Rat DRG Neurons	Lonza	R-DRG-505
PNGM™ BulletKit™	Lonza	CC-4461
NGF- β Nerve Growth Factor Beta	Sigma-Aldrich	N2513
Uridine	Sigma-Aldrich	U3003-5G
5-Fluoro-2'-Deoxyuridine	Sigma-Aldrich	F0503-100MG
50% Polyethylenimine Solution (PEI)	Sigma-Aldrich	P3143
Dulbecco's PBS without Ca ²⁺ /Mg ²⁺	Thermo Fisher	14040
Laminin	Sigma-Aldrich	L2020
Pipettes and Pipettors	Various	
15 mL and 50 mL Centrifuge Tubes	Various	

Equipment

Item	Vendor	Catalog #
Maestro MEA System	Axion BioSystems	
Axion Integrated Studio (AxIS)	Axion BioSystems	
37°C Water Bath	Various	
Cell Culture Incubator	Various	
Hemocytometer or Cell Counter	Various	
Biological Safety Cabinet	Various	
Tabletop Centrifuge	Various	
Phase Contrast Microscope	Various	
Liquid Nitrogen Storage	Various	