

MAESTRO APEX™

Automation for every facet of Maestro™ experiments

Introducing Maestro APEX. The all-in-one, plug-and-play automation solution for all aspects of MEA plate preparation, maintenance, and experiment execution. Seamlessly incorporating Hamilton robotics and a 44-plate capacity automated incubator with the Maestro™ platform and AxIS™ software, this integrated system provides a sterile environment for all steps leading up to and including running your Maestro experiment.

Have a lot of MEA plates to prepare?

Let APEX coat, wash, and spot cells onto your MEA plates. Load reagents and walk away knowing your plates will be prepared consistently and accurately.

Need to keep track of your plates?

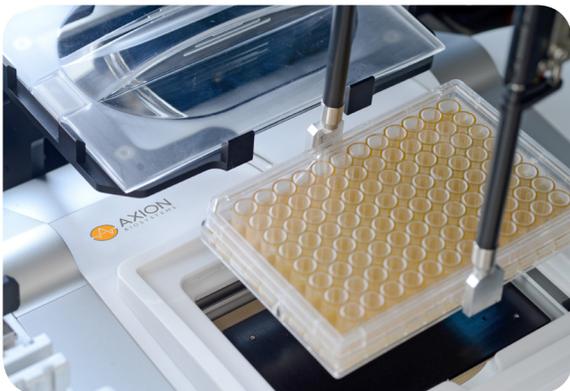
APEX includes two barcode readers, one embedded in the incubator and the other on the robotic arm.

Tired of endless hours in the hood maintaining your cell cultures?

The automated incubator and custom robotics protocols can perform media exchange at any time - including weekends.

Want to work off-line?

The Maestro can be removed for benchtop operation and later reinstalled onto the robotic plate deck with ease.



APEX PERFORMS

MEA Plate Preparation

- Matrix deposition
- Plate washing
- Cell spotting
- Media addition

MEA Plate Maintenance

- Media exchange

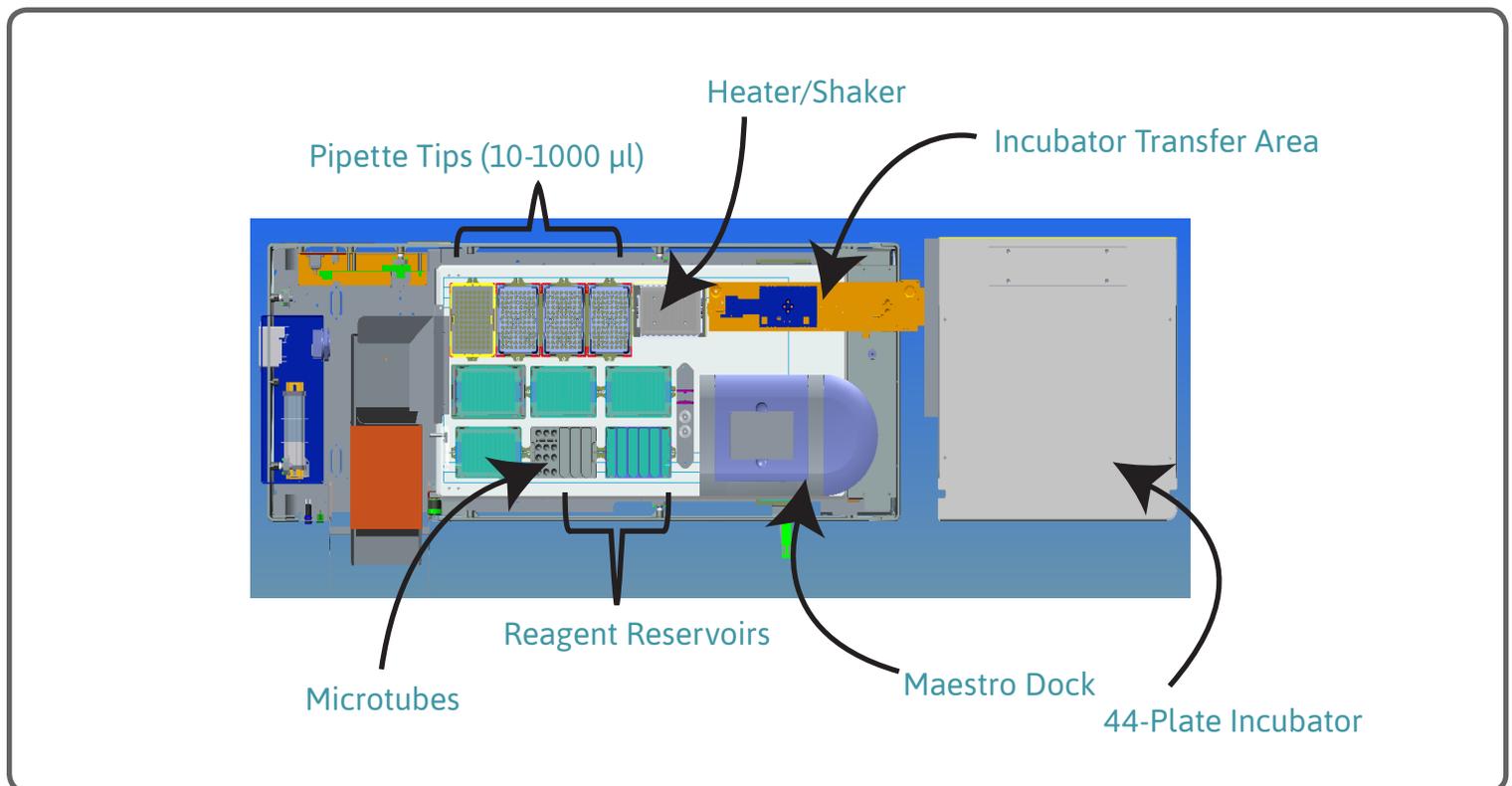
Maestro Experiments

- MEA plate docking
- Compound delivery
- CO₂ mixing and delivery



ROBOTIC DECK LAYOUT

(L x W x H = 74" X 28" X 52")



Feature

Specification

Sterile Plate Deck	HEPA filter and UV light package included
Integrated Incubator	Liconic, 44-plate capacity, automated plate handling; temperature control 33-70° C; 95% max humidity
4-Channel Pipettor	Fits 10-1,000 µl tips, transfer 1 µl-1,000 µl, volume verification at installation; exchangeable grippers for movement of plates
Reagent Reservoirs	50-60 ml capacity; standard deckware holds up to 9
Pipette Tip Racks	Holds 10 µl, 50 µl, 300 µl or 1,000 µl tips; houses up to 10 racks
Biohazard Waste Disposal	Solid waste disposal port and container for tips; 2-gallon liquid waste container
Barcode Readers	One each on the plate deck gantry and inside the incubator
ECmini™	Updated design providing responsive CO ₂ delivery; accommodates top-access to the MEA plate for the pipettor
Software	Complete platform operational control (robotics, incubator, AxIS) with Hamilton scheduling software; training included
Warranty	1-year full warranty

