

VIROLOGY STUDIES

SIMPLER, FASTER ASSAYS FOR INFECTIOUS DISEASES

Tired of waiting weeks to see results from plaque assays?

Annoyed by labor intensive and low-throughput endpoint assays methods?

Would you like to reduce the exposure to viral pathogens for your lab personnel?

Want to see reliable data faster and easier?

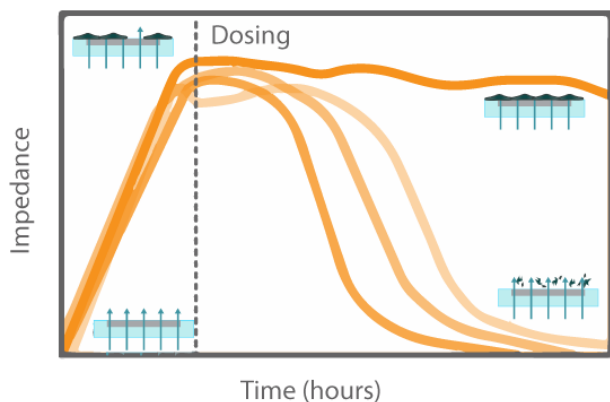
INTRODUCING THE MAESTRO Z...

Virus-induced cytopathic effects (CPEs) are usually associated with changes in cell morphology, strength of cell-cell contacts or cell-substrate attachment, and cell viability. Axion BioSystems' Maestro Z impedance platform can quantify all these dynamic cellular changes and provide real-time data analysis, revealing the kinetics of cell and virus interactions for improved mechanistic understanding.



APPLICATIONS

- Virus Titers
- Neutralization Assays
- Anti-viral Drugs
- Viricidal Efficacy
- Vaccine Development
- Oncolytic Virus Development



ADVANTAGES FOR VIROLOGY STUDIES

- **Label-free, non-invasive** impedance recording for long term monitoring.
- **It's easy!** Just culture your cells in the CytoView-Z plate add your treatment/therapy, and use the Maestro Z to monitor the killing kinetics in real time.
- Axion's 96-well CytoView-Z plates facilitate **high throughput** data collection and viewing of your cells under the microscope.
- AxIS-Z software eliminates labor intensive manual plaque counting by providing **real-time and automated data analysis**.
- **Environmental controls** provide a stable environment for both acute and long-term studies.



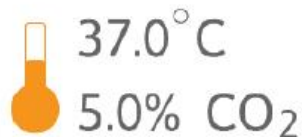
MAESTRO Z



96
ELECTRODES



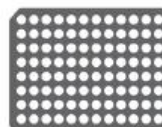
0.5 TB storage space



37.0° C
5.0% CO₂
Integrated Environment Chamber



Turn-key data analysis software



96- well assay plate



Viewing window for cell visualization