

AtlaZ

Virology research

Advance your virology studies

Using sophisticated impedance spectroscopy methodology, AtlaZ will allow you to accelerate your therapy and vaccine development by allowing monitoring of viral cytopathic effects (CPE) in real time, label-free and in high throughput format.

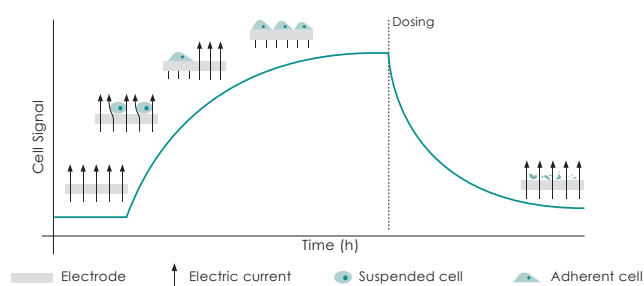
In order to gain a deeper mechanistic understanding of viral effects on cellular morphology, inter-cellular contacts or cell viability, real-time and continuous monitoring is necessary to access kinetic and phenotypic data.

Impedance assay applications:

- oncolytic virus development
- antiviral drug development
- vaccine development
- virucidal efficacy
- monitoring of viral cytopathic effects

Key benefits for virology research:

- **Real-time, label-free, long-term** impedance recordings (Cell Signal).
- **High throughput** with 6 x 96 well plates.
- **Electrical Impedance Spectroscopy**
- **Pre-set software modules** for easy calculation of cytolysis and Kill Time 50 values.
- **Automated data analysis** with access to raw data.



Contact us today



AtlaZ

Next level of live-cell analytics

AtlaZ Control Unit

- controls AtlaZ recording unit
- equipped with status display reflecting recording status in each of the 6 plates
- connects the recording unit with the laptop computer:
 - easy to use software with access to raw data
 - automated graphing of results
 - 21 CFR Part 11 compliance in GLP/GMP labs (coming soon)

AtlaZ Recording Unit

- 576 amplifier channels
- up to 6 x 96 plates simultaneously or independently
- real-time cell recording and analysis
- label-free experiments
- electrical impedance spectroscopy
- physiological conditions
- transparent 96-well plates
- barcode reader



Applications

- Immuno-oncology / CAR T
- Cytotoxicity
- Virology
- Cell characterization / QC
- GPCR / Receptor signaling
- Barrier function (TEER)
- Cell adhesion
- Wound healing