

AtlaZ

Quantitative live-cell analytics with GxP data compliance

Data integrity for impedance assays in GxP environment

For researchers seeking precise, real-time insights into cellular behavior, **AtlaZ** sets a new benchmark in live-cell analytics. By leveraging electrical impedance spectroscopy, AtlaZ provides a comprehensive, label-free approach to studying cell adhesion, proliferation, and cytotoxicity; GPCR signaling and cell morphology and barrier function - all in real time.

AtlaZ isn't just powerful - it's intuitive. **AtlaZ Control** software is designed for simplicity without compromising functionality. Its standardized measurement modes include pre-designed recording and analysis settings, ensuring a seamless workflow from experiment setup to data analysis.

All this information is presented in a single screen, allowing for efficient real-time decision-making and streamlined analysis. **FDA 21 CFR Part 11** compliance is available with a dedicated module.

Why choose AtlaZ system?

- **Real-time, label-free, long-term** impedance recordings
- **High throughput** with 6 x 96 well plates
- **Electrical Impedance Spectroscopy**
- **Continuous cell monitoring** to track cellular changes across a **broad frequency range**
- **Pre-set software modes** for easy calculation of cytolysis and Kill Time 50 values; IC_{50}/EC_{50} calculations to determine **drug potency**
- **Automated data analysis** with access to raw data
- **FDA 21 CFR Part 11 compliance** in GxP labs



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
 - FDA 21 CFR Part 11 compliance in GLP/GMP labs

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



AtlaZ GxP Software Module

- FDA 21 CFR Part 11 compliance
- system access and user tracking
- comprehensive audit trail
- tracked information includes:
 - unique data Identifier
 - measurement status
 - date & time stamps
 - user identification
- data integrity and protection:
 - immutable data handling.
 - audit-protected modifications
 - prevention of data tampering