

AtlaZ GPCR research

In depth GPCR studies

The AtlaZ system offers a highly automated approach capable of studying the pharmacology of various GPCR signaling pathways using a single platform. It requires no labeled reagents and provides the cell response to receptor stimulation in real-time.

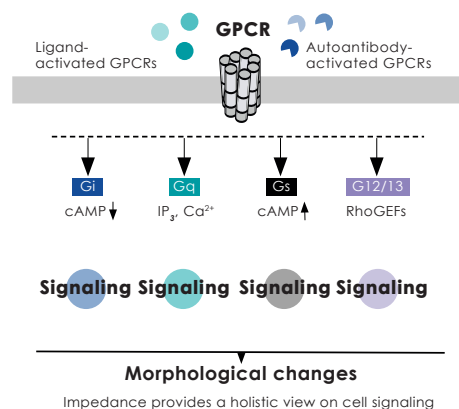
Impedance-based monitoring of GPCR stimulation is sensitive enough to work with endogenously expressed receptors in both primary cultures and stem cells.

Impedance assay applications:

- pharmacology of GPCRs
- monitoring of major G-protein-dependent pathways
- GPCR-mediated signal transduction effects
- endogenously expressed receptors
- detection of morphological changes of cells

Key features for GPCR research:

- **Real-time, label-free, long-term** impedance recordings (Cell Signal)
- **High throughput** with 6 x 96 well plates
- **Electrical Impedance Spectroscopy (EIS)**
- **Continuous** cell monitoring
- **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)

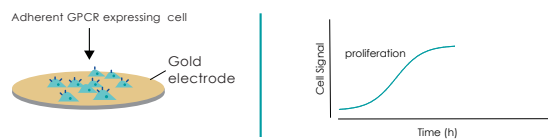
AtlaZ Recording Unit

- 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

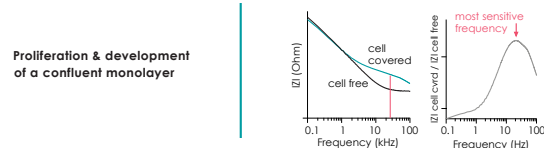


Workflow of a GPCR assay

1 Cell plating and proliferation



2 Find optimal EIS frequency



3 GPCR profiling

