

# AtlaZ

## Immuno-oncology research

### Kinetic effects, label-free and in real-time

Real-time impedance recordings provide insights in various cancer cell phenotypes, such as cell morphology, proliferation, lateral migration, as well as cytotoxicity by biologics or chemical entities over prolonged periods of time.

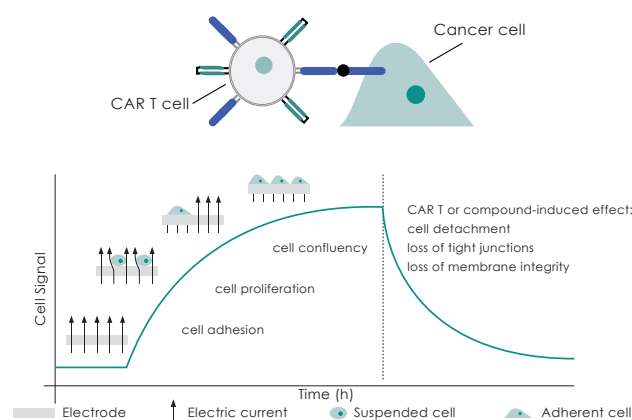
A crucial advantage over standard assays is the continuity of cell monitoring. Such measurements reveal the kinetics of cell behavior and allow an in-depth mechanistic understanding without the need for time- and labor-intensive endpoint assays. For immuno-oncology assays, this means easy in-depth insight into potency and kinetics of immune cell killing of cancer cells.

#### Impedance assay applications:

- Immuno-oncology assays
- CAR T cell potency assessment
- NK cell potency assessment
- Ab-dependent cellular cytotoxicity (ADCC)
- Liquid tumors cell assay

#### Key benefits for immunology 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 cytotoxicity 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