

Is your safety program up-to-date? Keep your pole position with Patchliner & CardioExcyte 96!



Patchliner.
High quality patch clamp screening.

Comprehensive cardiotoxicity screening

CardioExcyte 96 reveals compound effects on beating parameters recorded from intact cardiac networks.

Patchliner allows detailed investigations on ion channel level. In combination, these platforms provide unmatched information on a compound's safety profile.

CardioExcyte 96 & Patchliner offer you:

- Efficient and reliable cardiotoxicity profiling and screening using stem cell-derived cardiomyocytes
- Compatibility with FDA-, HESI- and Task Force Five-stipulated requirements
- Excellent compatibility with all commercially available, pluripotent stem cells



CardioExcyte 96.
Label-free impedance system for
beating cardiac networks.

CardioExcyte 96 & Patchliner

nanjion

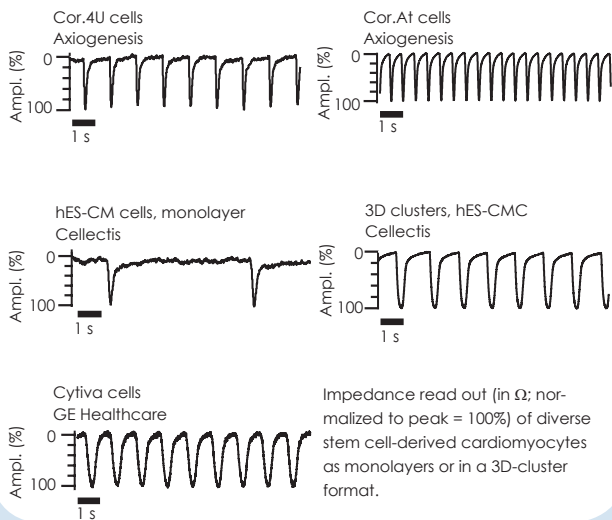
Fast and reliable cardiac safety screening

"Late compound withdrawal due to cardiac safety liability issues is costly and inefficient. Platforms are required supporting reliable compound testing in early development. The Patchliner has drastically improved the efficiency of our safety testing by applying a high degree of automation to otherwise manually performed low throughput patch clamp experiments with maintained data quality. The compatibility with stem cells is extraordinary, a vital feature regarding the upcoming need for the Comprehensive In vitro Proarrhythmia Assay (CIPA) recently discussed at the FDA/CSRC/

HESI Think Tank Meeting. The new paradigm aims at mechanistic understandings of ventricular arrhythmias by assessing multiple human cardiac currents and action potentials in stem cell-derived cardiomyocytes in combination with in silico modeling of ventricular myocyte physiology. The combination of cardiac network measurements followed by high quality patch clamp increase the possibility substantially, to quickly and correctly assess a compound's cardiotoxicity profile"

Dr. Liudmila Polonchuk, Principal Scientist, Non-Clinical Safety, F. Hoffmann La Roche

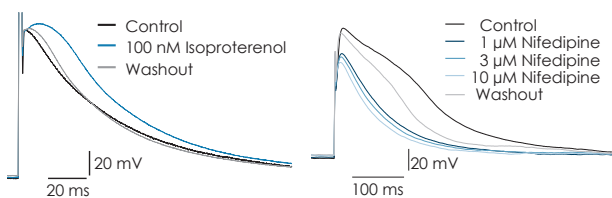
Beating cardiomyocyte networks



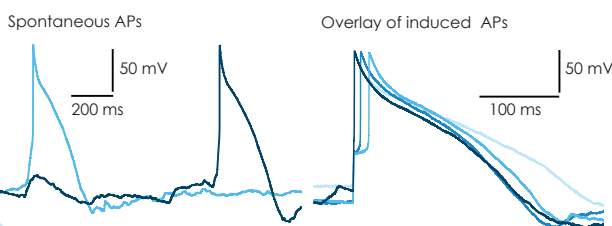
CardioExcyte 96:

- Non-invasive, label-free measurements of beating cardiomyocyte networks
- 96 recording wells in parallel with 1 ms time resolution
- Quick experiments or long-term compound effects on cardiotoxicity
- Real-time access to beating parameters
- Outstanding software for data analysis and export
- Cost efficient consumables - 96-well format

Pharmacology on action potentials



Spontaneous and induced action potentials



Patchliner:

- High quality patch clamp recordings of compound effects on individual ion channels
- Reliable and efficient compound screening on cardiac ion channels incl. hERG, Nav1.5 & L-type calcium
- Detailed investigations of compound effects on action potentials
- Recordings at physiological temperatures
- Cost-efficient consumables
- Powerful and efficient data analysis
- Up to 500 data points per day

CardioExcyte 96 & Patchliner. Pole position in cardiotoxicity screening.