

GREEN CHEMISTRY, RED FLAGS:

Multiparametric cardiotoxicity screening of phytochemicals using hiPSC-CMs-MEA Assay

● Monday, 27th of November 2023

● Laura-Sophie Frommelt



Importance of pre-clinical comprehensive cardiac safety studies

Cardiotoxicity:

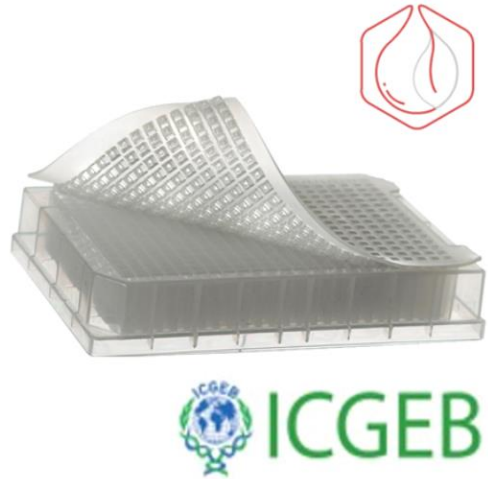
- is one of the most common forms of **drug-induced toxicity**
- Resulted in numerous costly **withdrawals of drugs**

Preclinical cardiac safety studies:

- Increase **drug safety**
- Reduce the **cost** of drug development

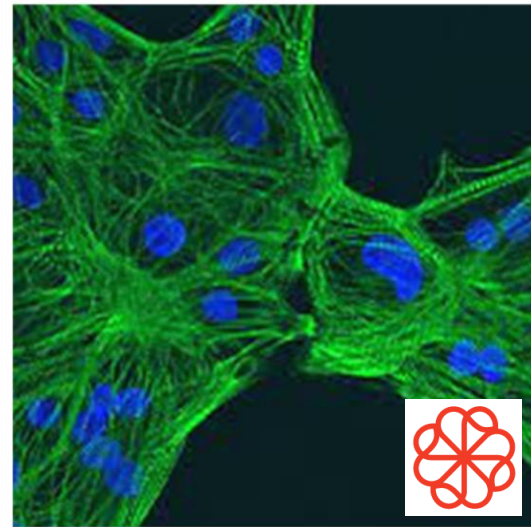
Aim: Assess the effectiveness of the **hiPSC-CMs-MEA assay** in detecting acute and chronic cardiotoxic effects of phytochemicals

Multiparametric drug assessment using hiPSC-CMs-MEA Assay



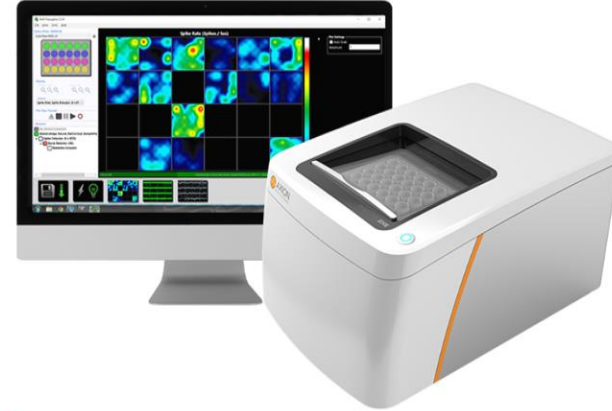
Prestwick Phytochemical Library:

- 320 purified well-known and novel compounds
- mostly from plants



Ncyte Cardiomyocytes (Ncardia)

- Human iPSC-derived ventricular-like CMs
- Large-scale production
- Slow + uniform beating rate

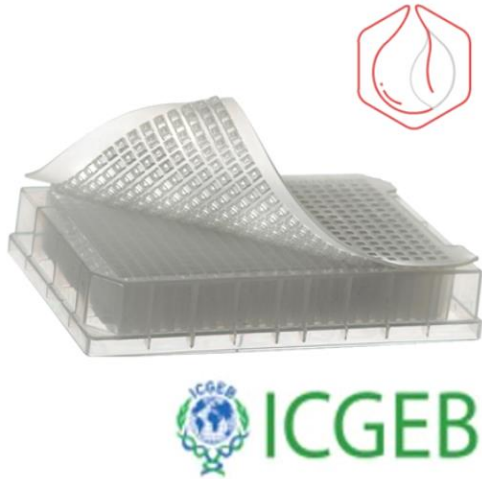


 **Maestro Edge MEA system**
Axion Biosystems

Maestro Edge MEA system:

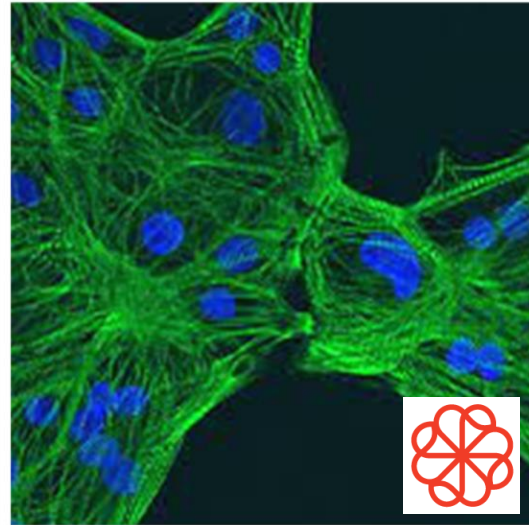
- CiPA-study (Comprehensive *in vitro* Proarrhythmia Assay)
- Heartbeat in a dish
- Various cardiac endpoints: FP-signals, LEAP, contractility signals and cell viability

Multiparametric drug assessment using hiPSC-CMs-MEA Assay



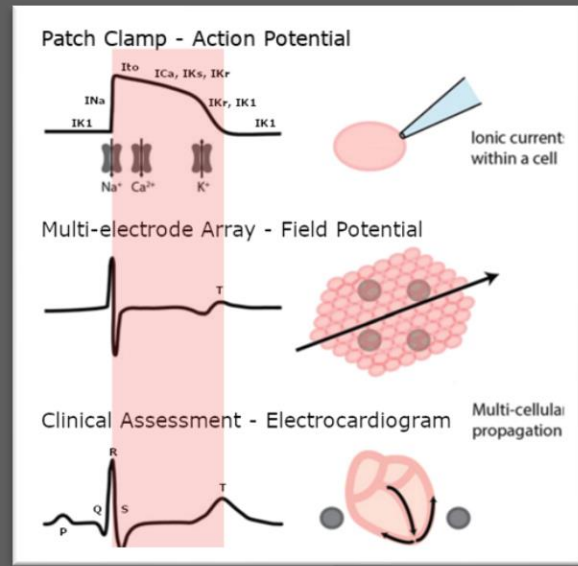
Prestwick Phytochemical Library:

- 320 purified well-known and novel compounds
- mostly from plants



Ncyte Cardiomyocytes (Ncardia)

- Human iPSC-derived ventricular-like CMs
- Large-scale production
- Slow + uniform beating rate

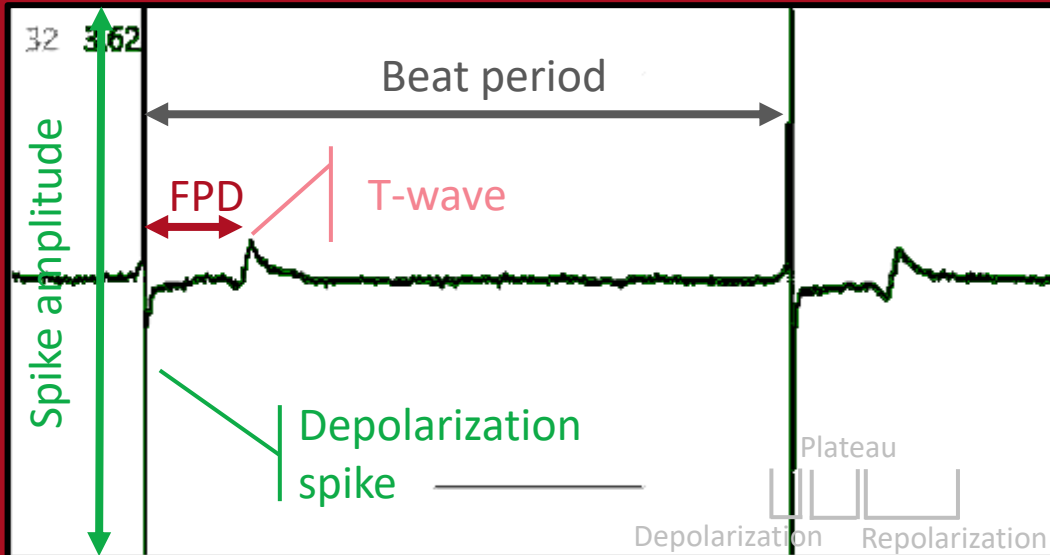


Maestro Edge MEA system:

- CiPA-study (Comprehensive *in vitro* Proarrhythmia Assay)
- Heartbeat in a dish
- Various cardiac endpoints: FP-signals, contractility signals and cell viability

Cardiac endpoint parameters

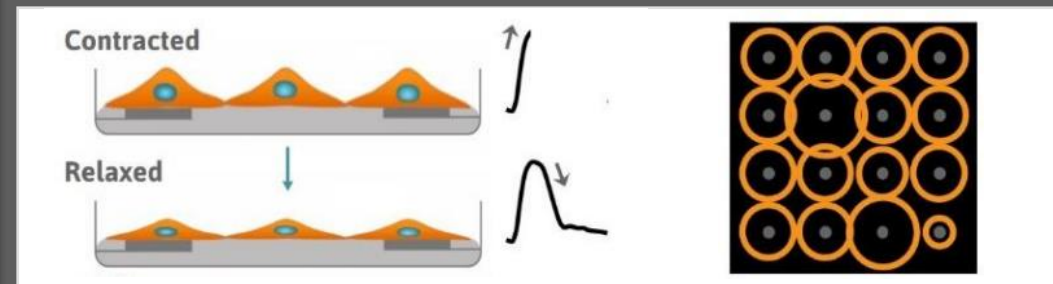
Field potential-based



- Beat period mean
- Beat period irregularities
- Field potential duration (+corrected)
- Spike amplitude mean
- Arrhythmic events
- Local extracellular action potential (LEAP)

Impedance-based

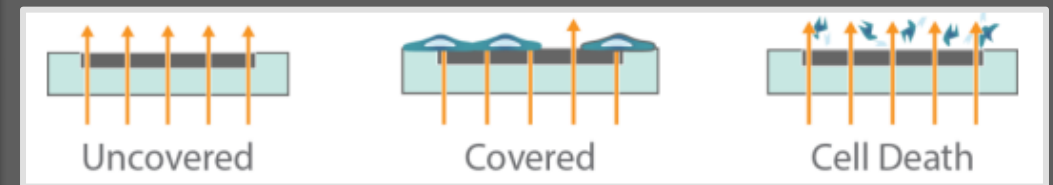
Contractility (3.125kHz)



Beat amplitude

Excitation-contraction coupling

Viability (41.5kHz)



Weighted mean resistance

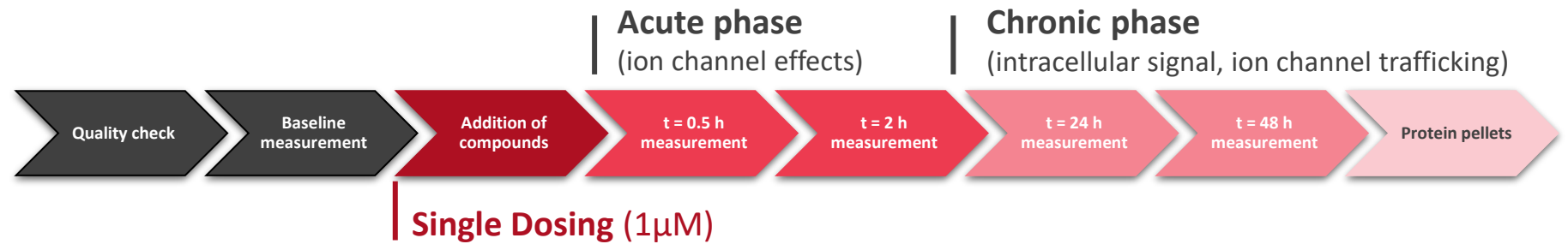
Average resistance

Experimental design of the library screening

Plate design

	1	2	3	4	5	6
A	Blue	Blue	Blue	Light Blue	Dark Red	Dark Red
B	Blue	Blue	Blue	Light Blue	Dark Red	Dark Red
C	Red	Red	Yellow	Yellow	Yellow	DMSO
D	Red	Red	Yellow	Yellow	Yellow	E-4031

Timeline



Positive control E-4031 (500nM)



Measurements:

- **Field Potential** (+paced)
- **Viability**
- **Contractility** (+paced)
- LEAP (2+24h)

Data analysis

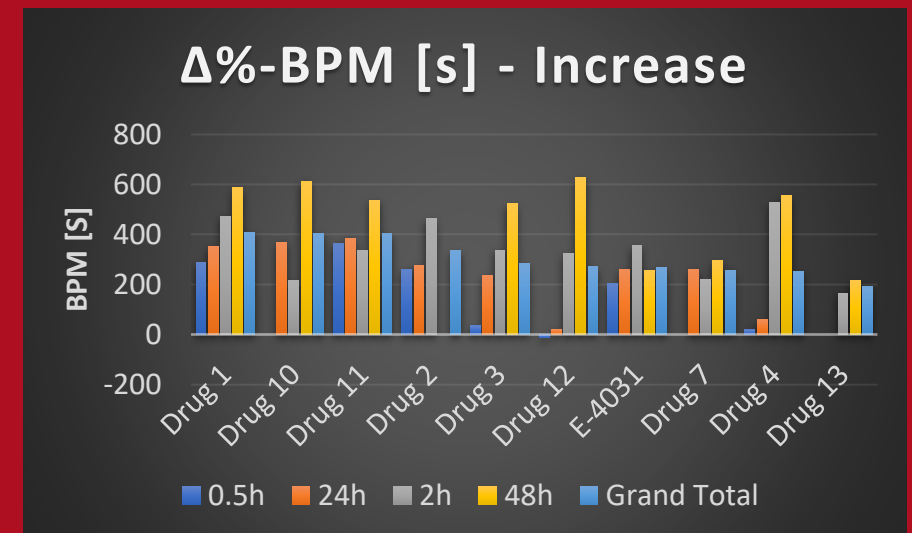
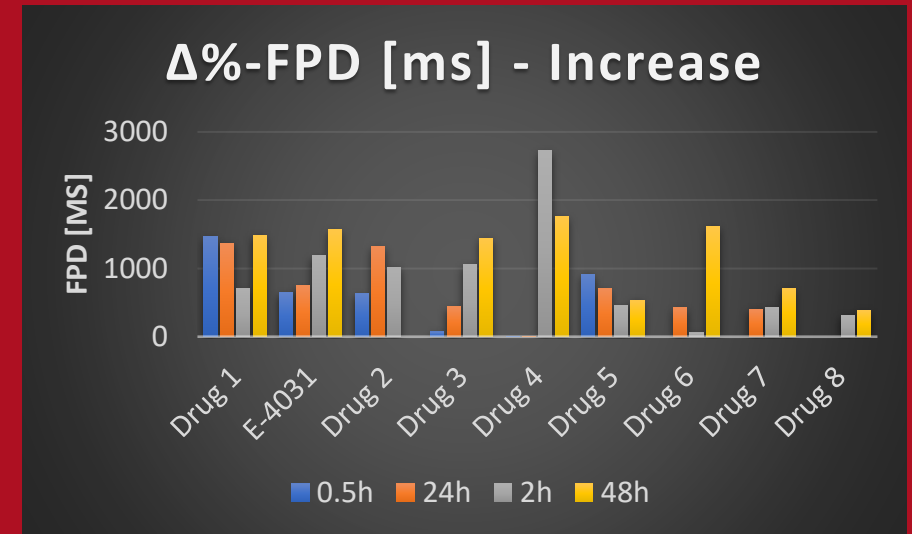
- **Normalized comparison:** %-change between baseline and dosed conditions
- **Subtraction** of the vehicle control (DMSO)
- **Min-Max Scalar normalization** \rightarrow **ToxPiTM Ranking**

Validation – K⁺- channel blockers (K_v11.1 - hERG)

E-4031 dihydrochloride (positive control)

Expectations

- Prolongation of Field Potential Duration (FPD)
- Increase in Beat Period Mean (BPM)
- Arrhythmic events

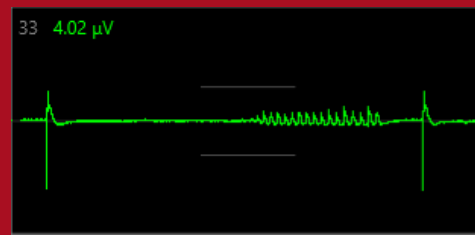
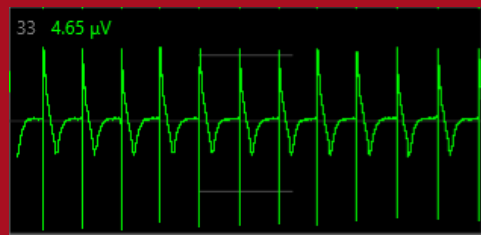


Validation – K⁺- channel blockers (K_v11.1 - hERG)

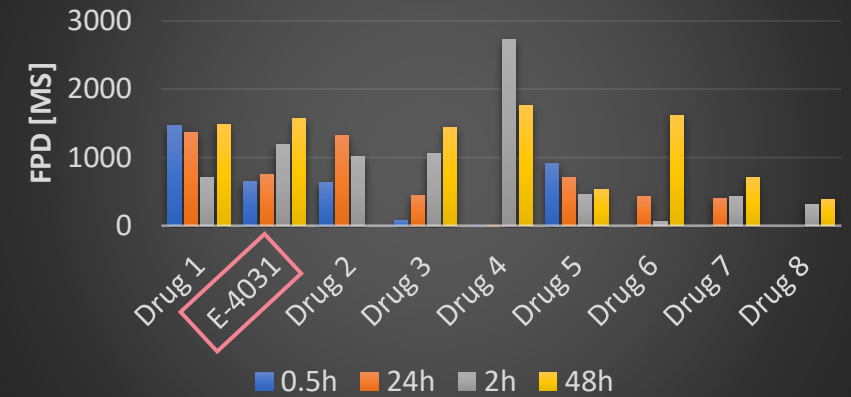
E-4031 dihydrochloride (positive control)

Results:

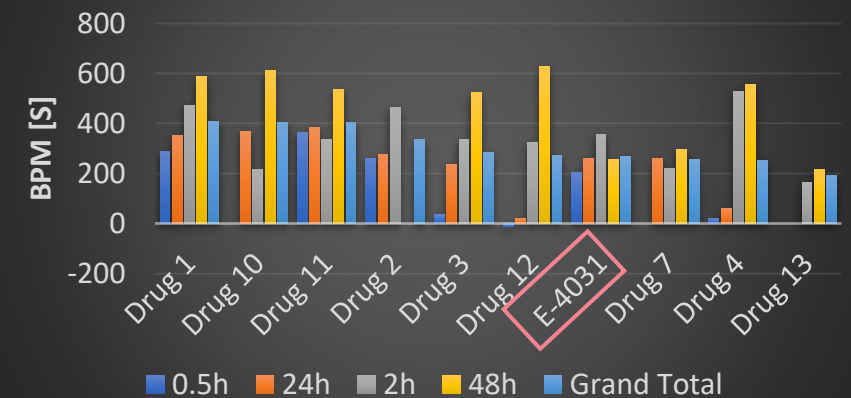
- Prolongation of Field Potential Duration (FPD): **2nd place**
- Increase in Beat Period Mean (BPM): **6th place**
- Arrhythmic events: **1st place (20 events)**



Δ%-FPD [ms] - Increase



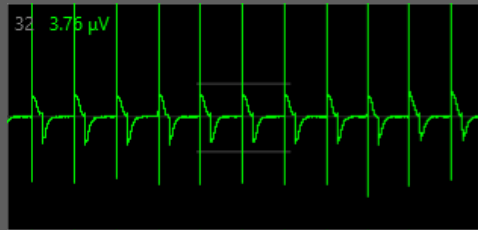
Δ%-BPM [s] - Increase



Validation – K⁺- channel blockers

Drug 1: Indole - Cardiovascular active compound, Vasodilator

- Arrhythmic events: 3rd place (17 events)



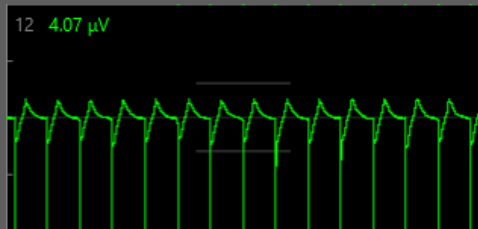
Baseline



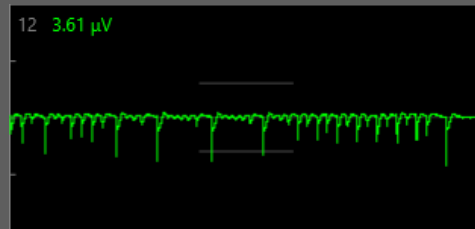
24h

Drug 4: Alkaloid - Cardiovascular active compound, Antihypertensive

- Arrhythmic events: 11th place (12 events)

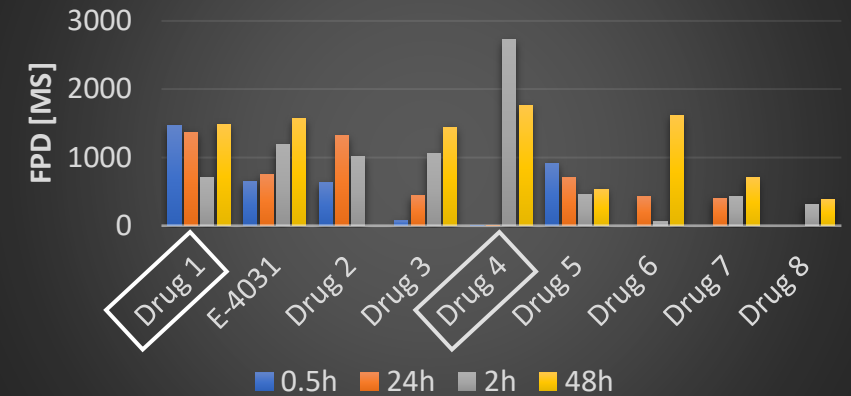


Baseline

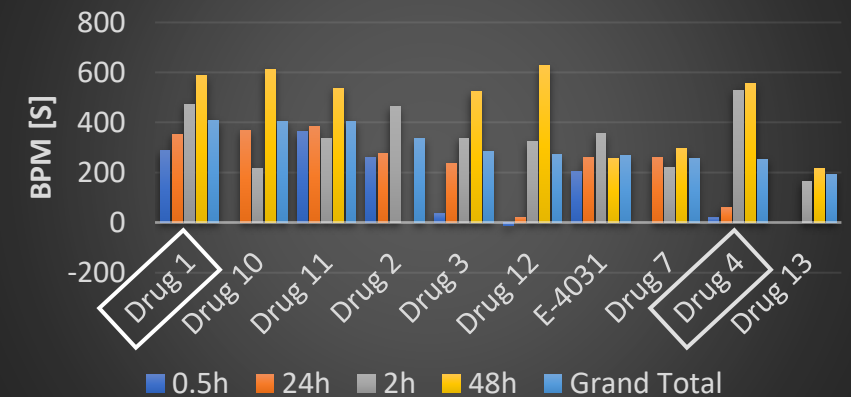


24h

Δ%-FPD [ms] - Increase

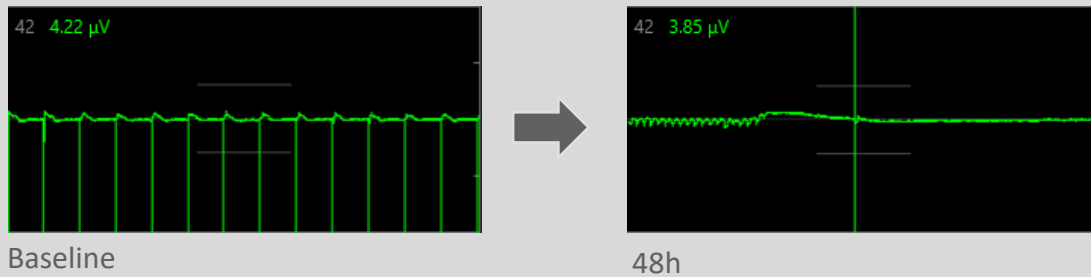


Δ%-BPM [s] - Increase

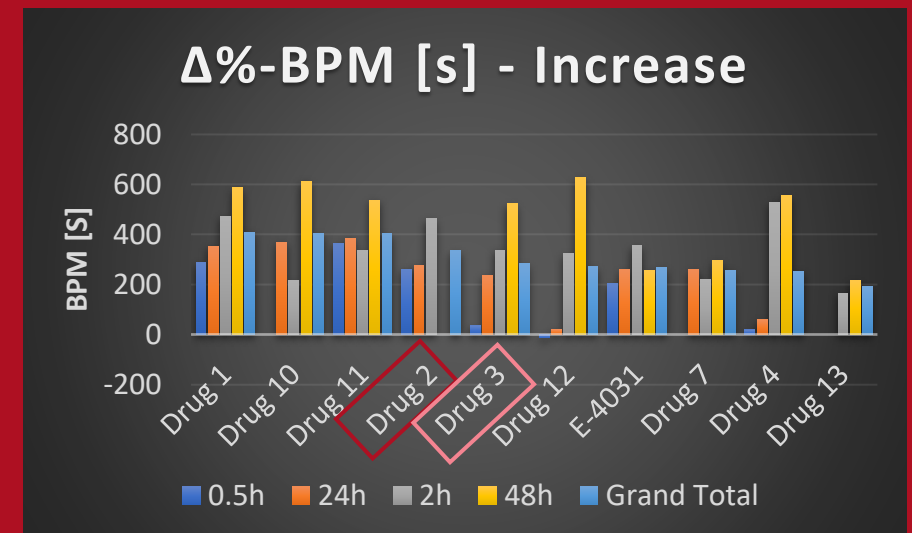
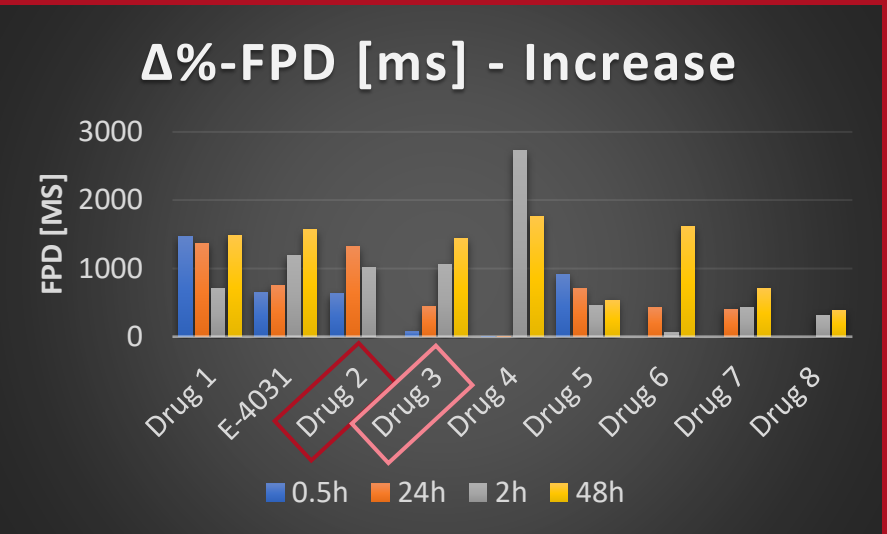
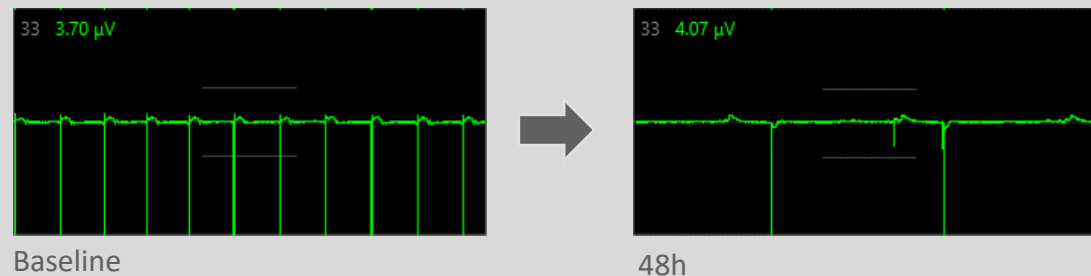


New Hits – possible K⁺- channel blockers

Drug 2: semisynthetic alkaloid – CNS active compound, Neuroprotectant
Arrhythmic events: 2nd place (19 events)

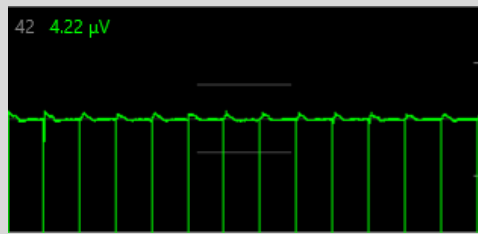


Drug 3: Isoquinoline, Infectiology, antibacterial effect
Arrhythmic events: 5th place (15 events)

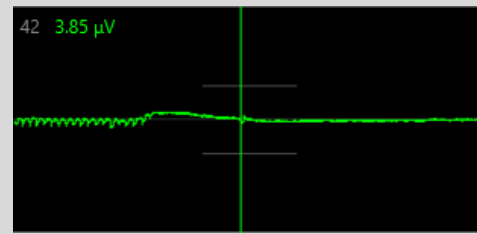


New Hits – possible K⁺- channel blockers

Drug 2: semisynthetic alkaloid – CNS active compound, Neuroprotectant
Arrhythmic events: 2nd place (19 events)



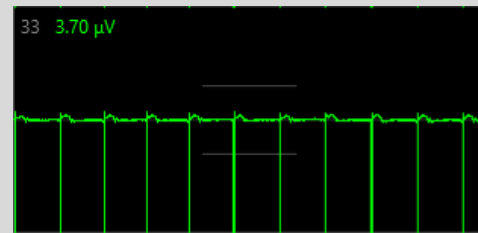
Baseline



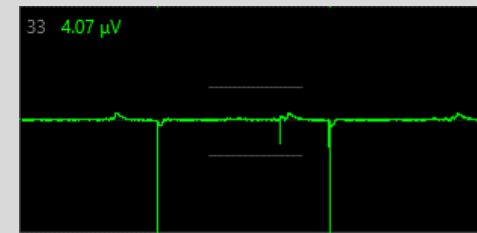
48h

Drug 3: Isoquinoline, Infectiology, antibacterial

Arrhythmic events: 5th place (15 events)



Baseline



48h

Δ%-FPD [ms] - Increase

3000

**NO reported link
with arrhythmia
so far**

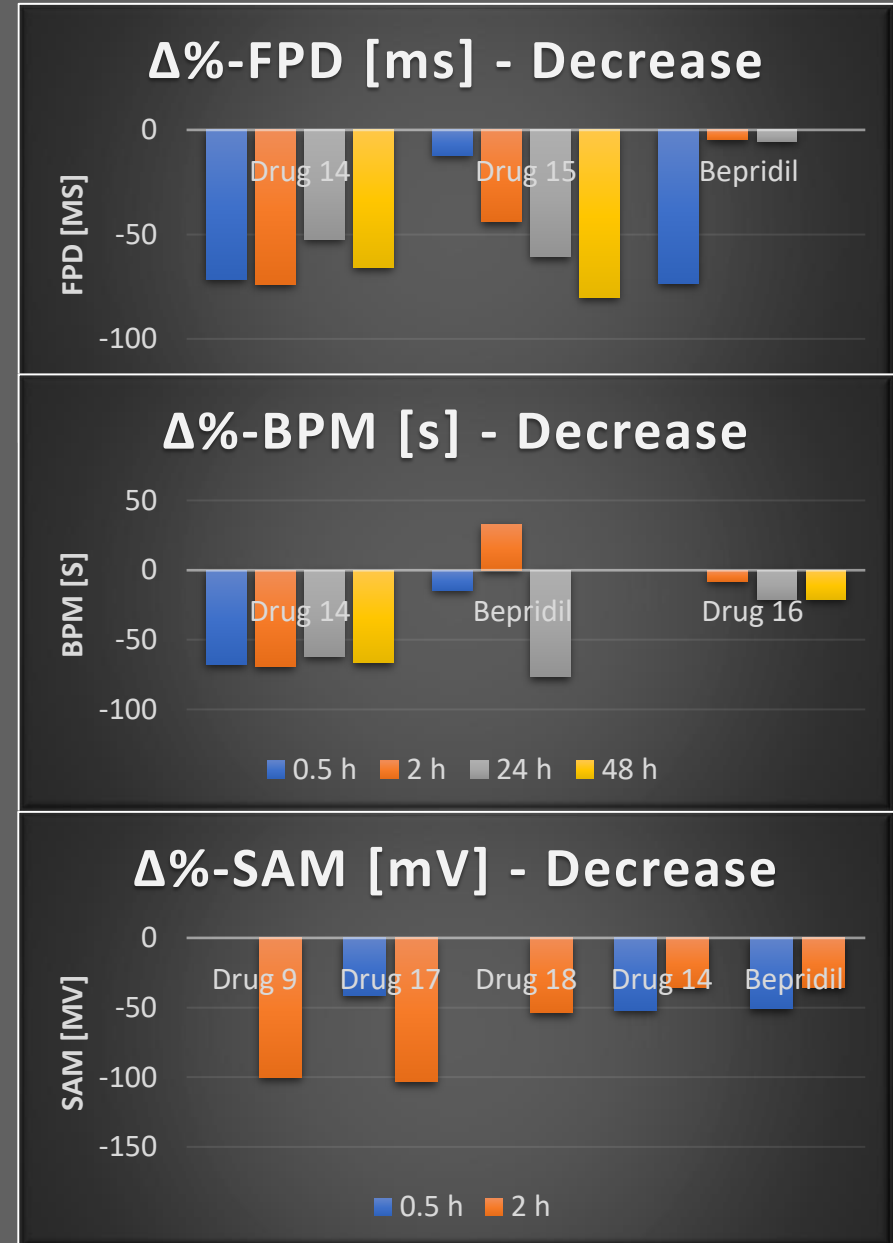


Validation – Na⁺/Ca²⁺-channel blockers

Bepridil (positive control) - Na⁺ and Ca²⁺-channel blocker, antiarrhythmic and antihypertensive

Expectations:

- Shortening of Field Potential Duration (FPD)
- Decrease in Beat Period Mean (BPM):
- Decrease in Spike Amplitude Mean (SAM)

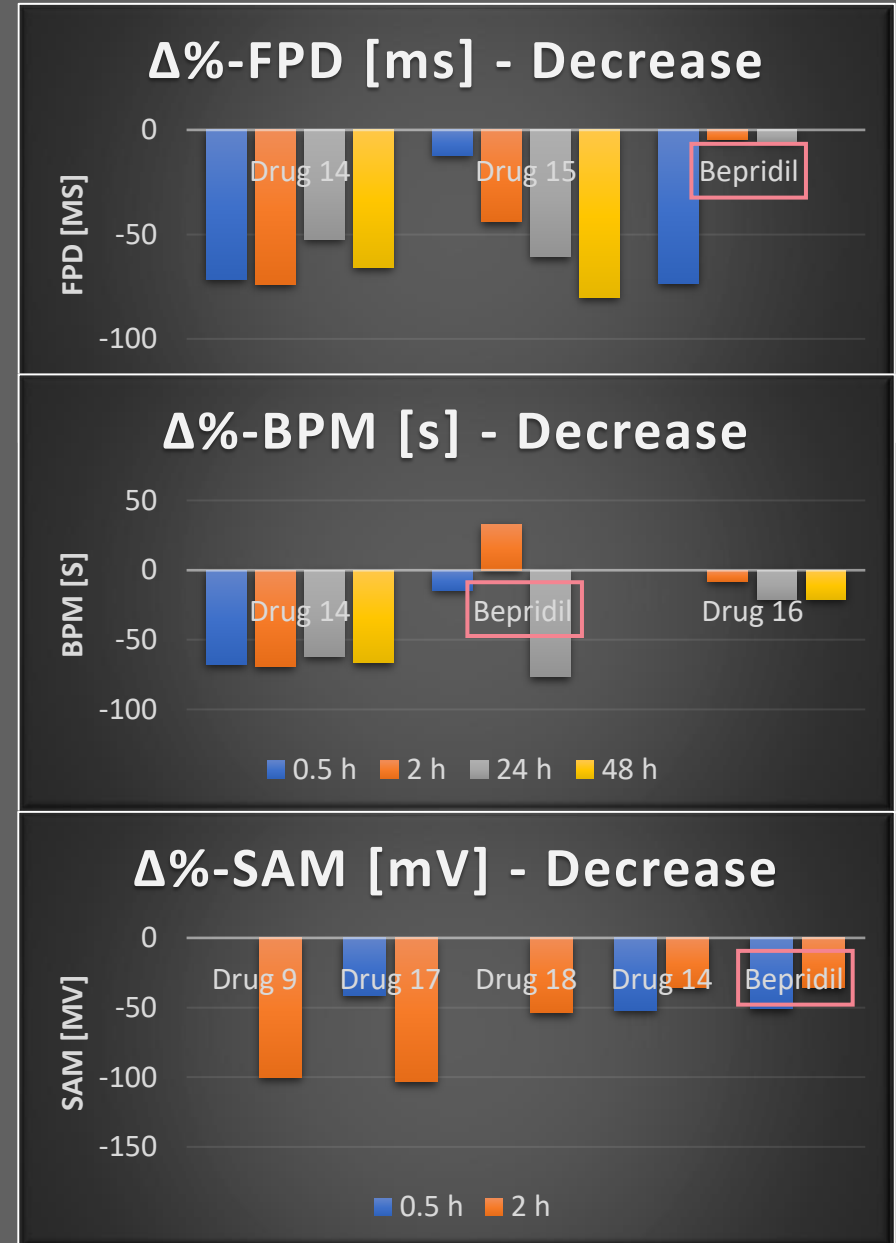


Validation – Na⁺/Ca²⁺-channel blockers

Bepridil (positive control) - Na⁺ and Ca²⁺-channel blocker, antiarrhythmic and antihypertensive

Results:

- Shortening of Field Potential Duration (FPD): **3rd place**
- Decrease in Beat Period Mean (BPM): **2nd place**
- Decrease in Spike Amplitude Mean (SAM): **5th place**

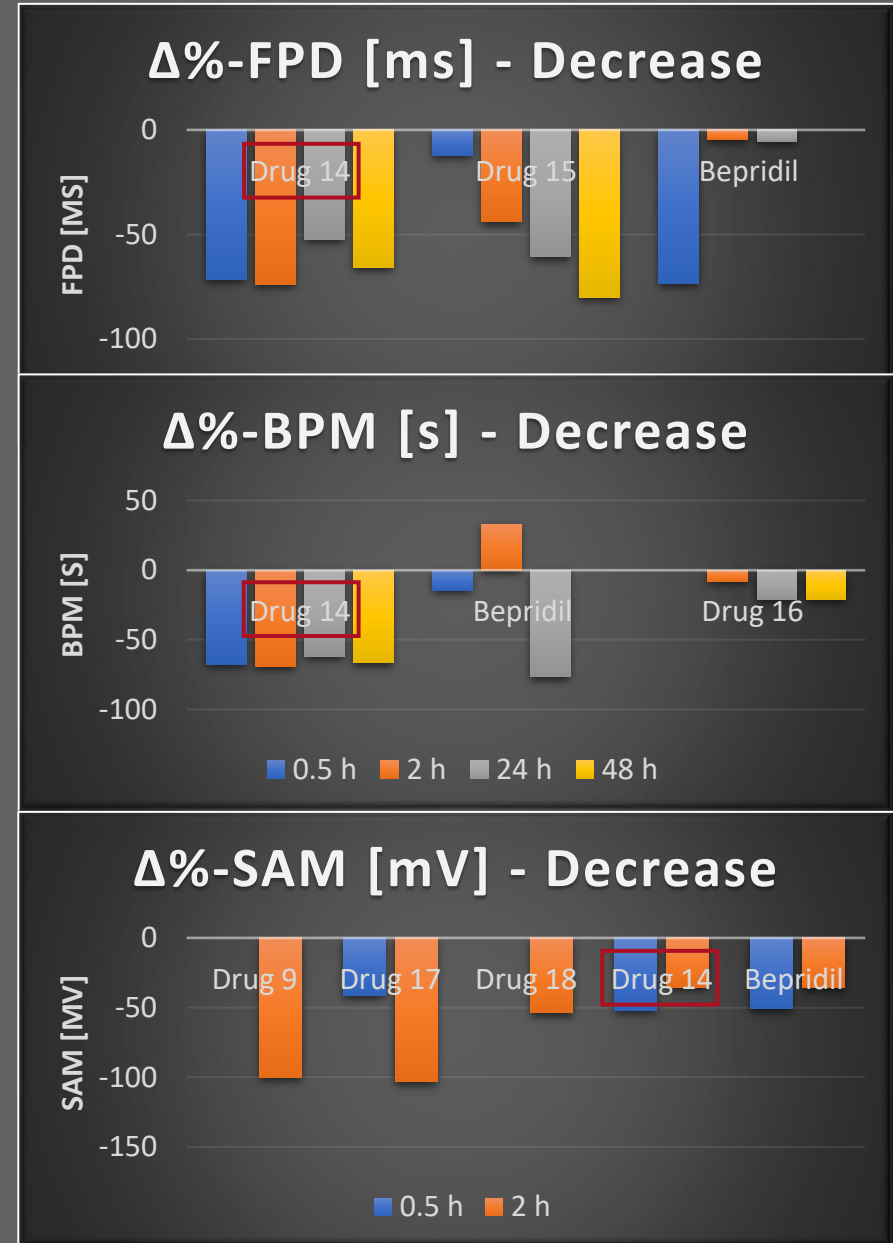


Validation – Na⁺/Ca²⁺-channel blockers

Drug 14: Alkaloid, Cardiovascular active compound, antihypertensive effect

- Shortening of Field Potential Duration (FPD): **1st place**
- Decrease in Beat Period Mean (BPM): **1st place**
- Decrease in Spike Amplitude Mean (SAM) : **4th place**

→ has been used in the treatment of hypertension but has largely been replaced by drugs with fewer adverse effects



Moving from single parameter to multiparameter analysis

Individual parameters:

Beat Amplitude: 69 compounds

Field Potential Duration: 13
compounds

Beat Period Mean: 17 compounds

Beat Period Irregularities: 23
compounds

Spike Amplitude Mean: 39
compounds

Severe arrhythmic effects: 27
compounds

Viability: 14 compounds

Multiparametric analysis aims to:

- quantify the **actual arrhythmogenic risk** rather than sole focus on a single parameter
- comprehensive characterization of cardiotoxicity
- early and informed decision:

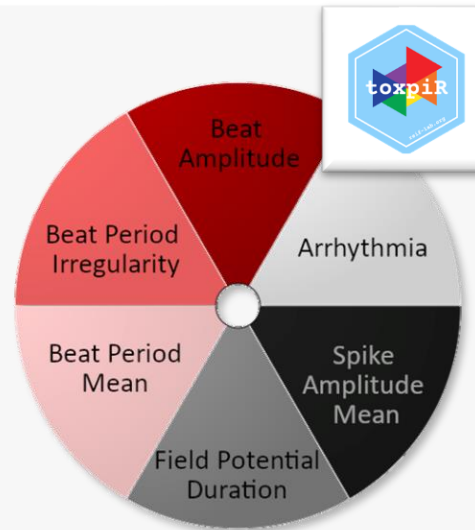
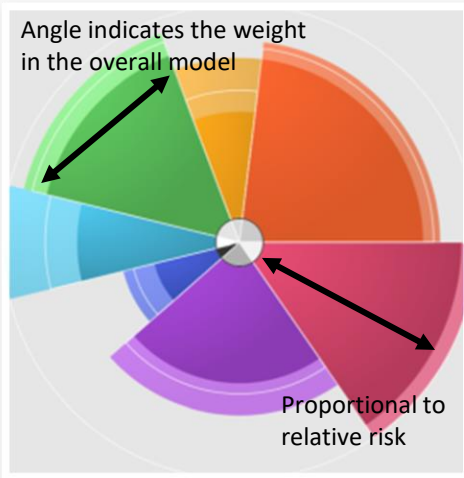


or



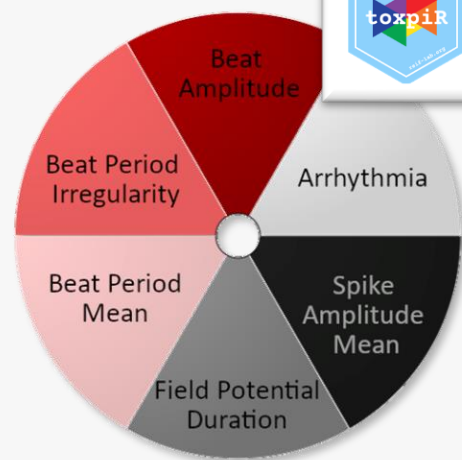
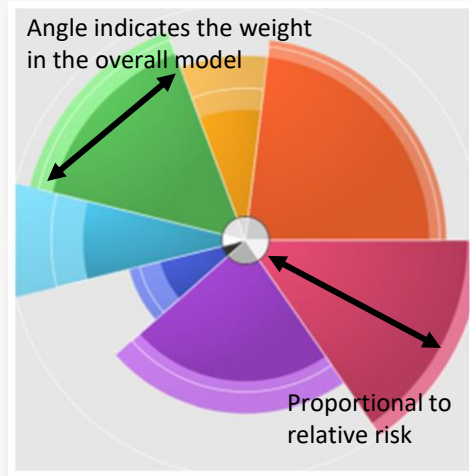
Moving from single parameter to multiparameter analysis

Beat Amplitude: 69 compounds
Field Potential Duration: 13 compounds
Beat Period Mean: 17 compounds
Beat Period Irregularities: 23 compounds
Spike Amplitude Mean: 39 compounds
Severe arrhythmic effects: 27 compounds
Viability: 14 compounds

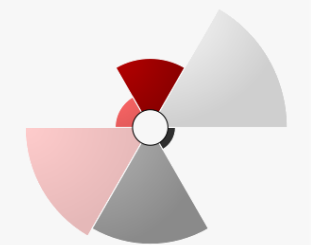
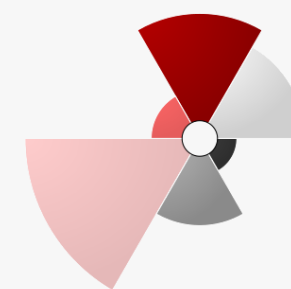
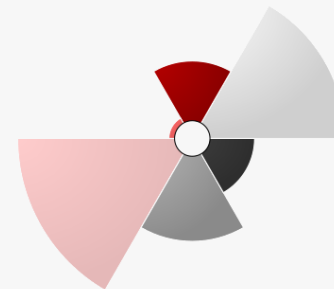
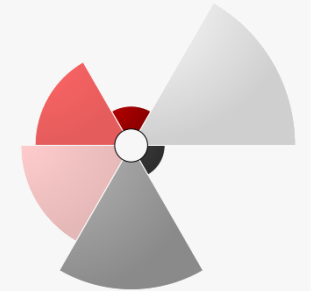
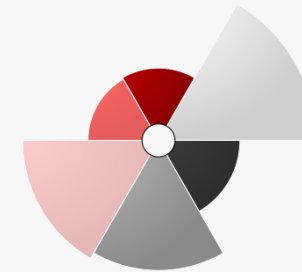
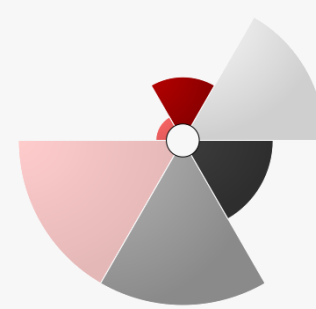


Moving from single parameter to multiparameter analysis

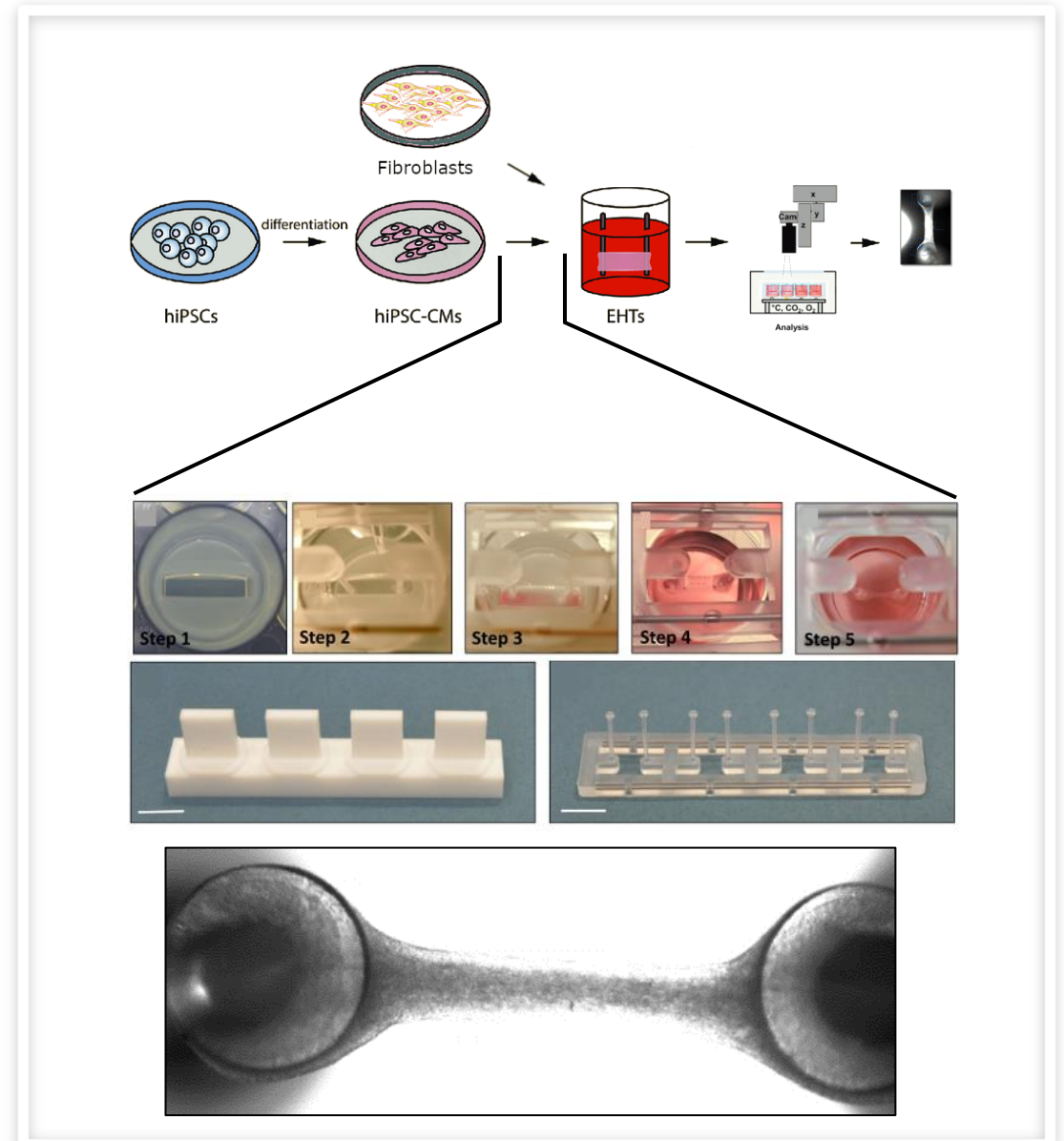
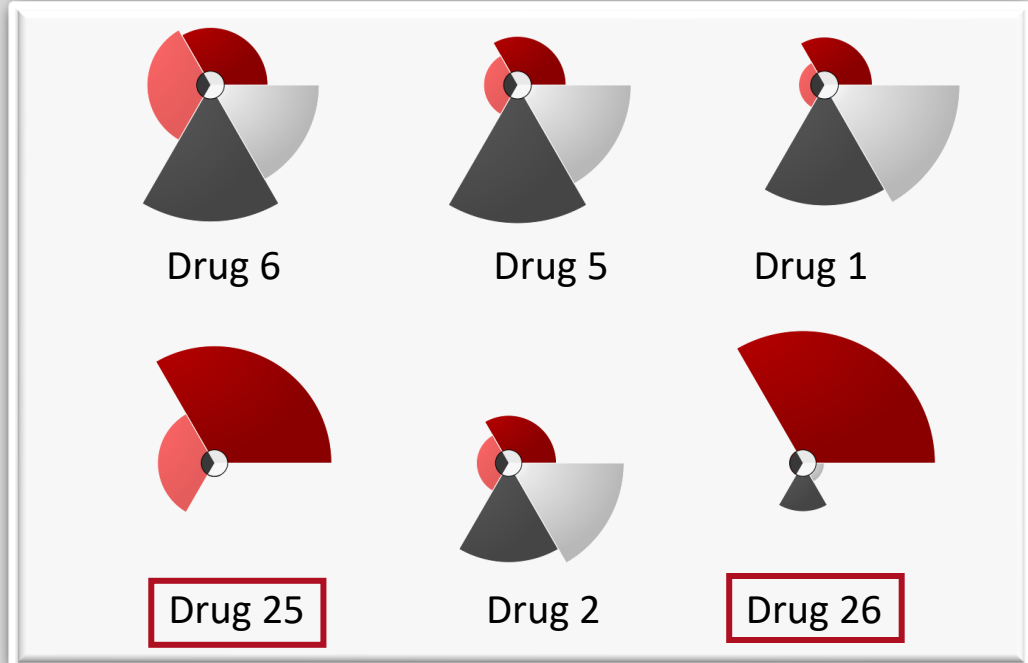
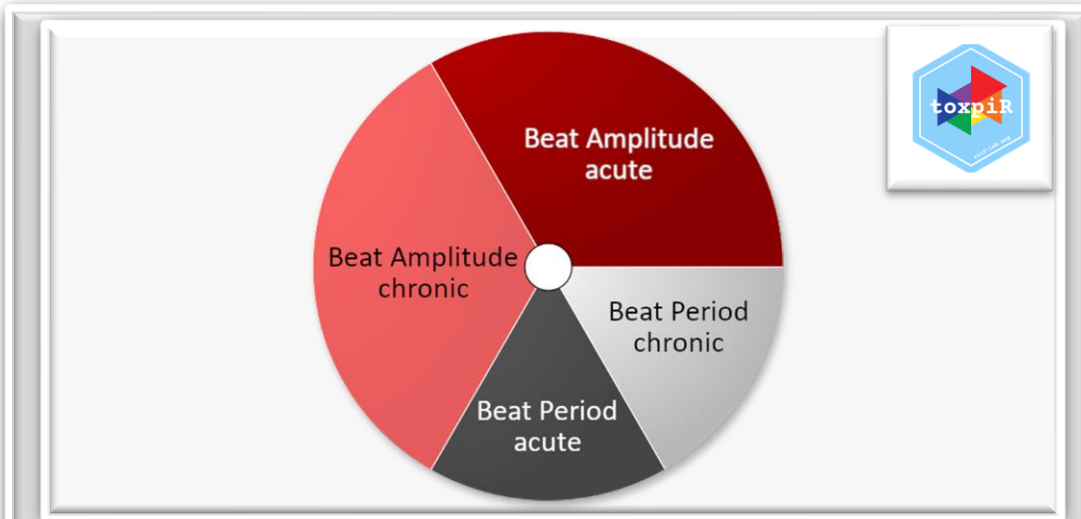
Beat Amplitude: 69 compounds
Field Potential Duration: 13 compounds
Beat Period Mean: 17 compounds
Beat Period Irregularities: 23 compounds
Spike Amplitude Mean: 39 compounds
Severe arrhythmic effects: 27 compounds
Viability: 14 compounds



Top six compounds:



Inotropic compounds



Summary

hiPSC-CMs-MEA Assay is a useful tool for preclinical assessment of cardiotoxic effects

- Positive controls (E-4031, Bepridile, etc.) are among the highest ranked compounds
- 81 compounds have a significant effect on 1 or more of the analyzed cardiac parameters
- Multiparameter analysis

eurac
research



Katarina Mackova



Serena Zacchigna



Giulio Ciucci



Alessandra Rossini



Chiara Volani

Acknowledgments

