

Next level toxicity screening: From single channel to overall cell behavior

nanijon

Conrad Weichbrodt¹, Mohamed Kreir¹, Matthias Beckler¹, Alison Obergrussberger¹, Ilka Rinke¹, Michael George¹,

Sonja Stoelze-Feix¹, Andrea Brüggemann¹, Niels Fertig¹

¹Nanon Technologies GmbH, Gabrielenstr. 9, 80636 Munich, Germany

Abstract

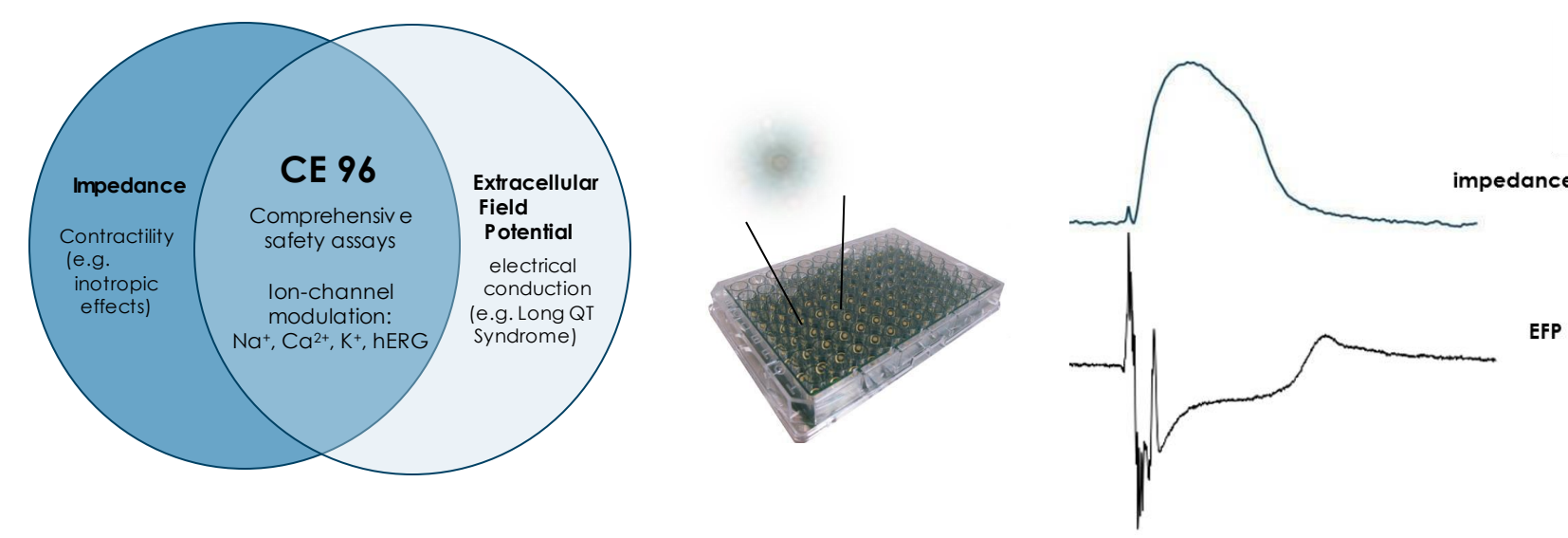
Many lead compounds fail in the late stages of the drug development process mainly by inflicting drug induced injury on liver, heart or other organs. Therefore, devices for detecting possible cell toxicity in early stages of the development process are highly demanded. Especially ion channels represent an important class of drug targets for in vitro pharmacological profiling.

High throughput screening (HTS) assays such as automated electrophysiological patch clamp and impedance based assays allow for the determination of drug effects on a whole cell level whereas artificial bilayers provide a robust environment for the assessment of single ion channel molecules.

We here present the CardioExcyte 96, a system providing a combination of Electric Impedance Spectroscopy (EIS) as well as Electric Field Potential (EFP) readout for a network of diverse cells like iPSC cardiomyocytes or hepatocyte-like cells which is exemplified by toxicity effects utilizing reference compounds such as Dofetilide (on iPSC cardiomyocytes) or Paracetamol (on hepatocyte-like cells).

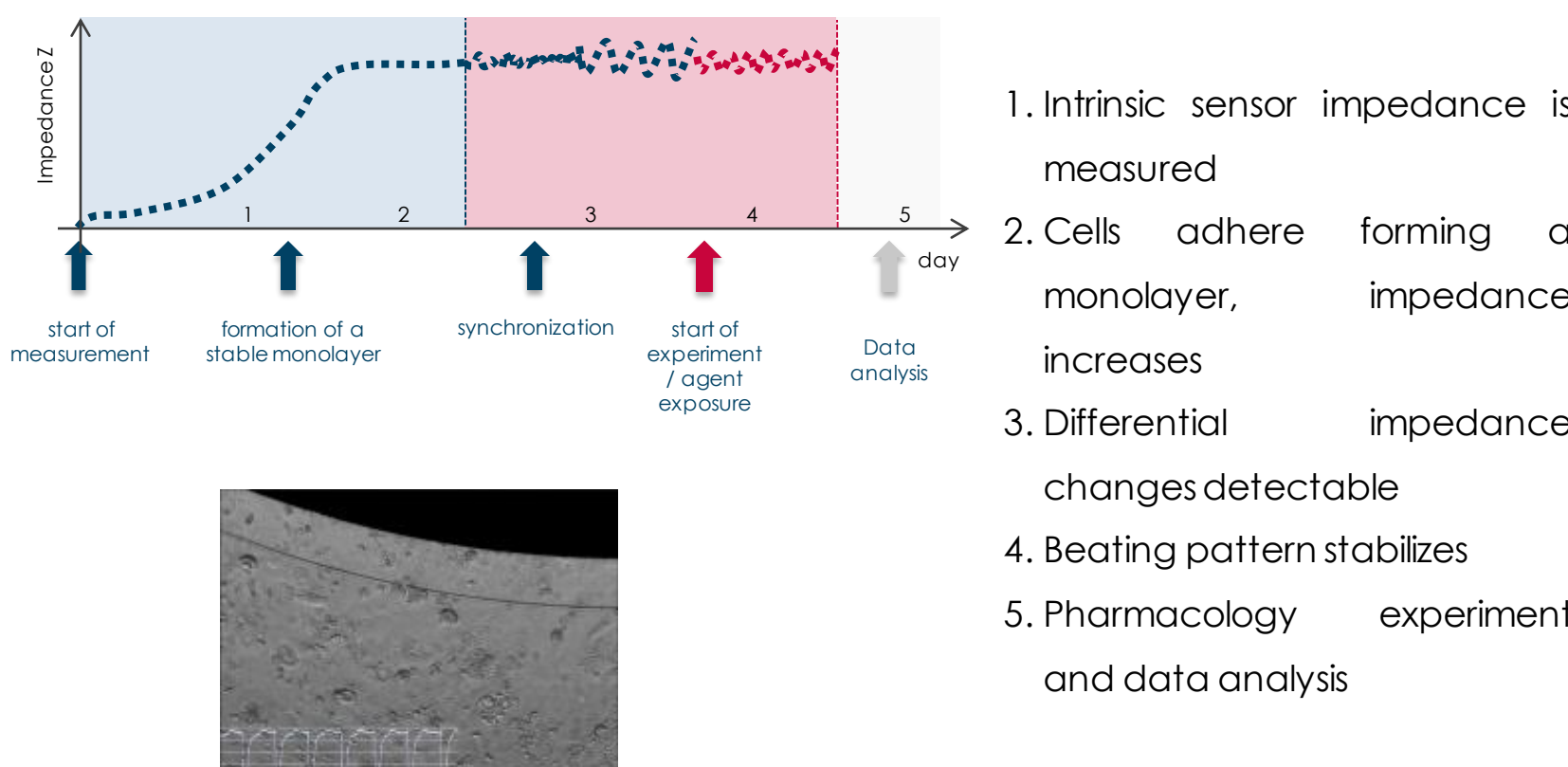
Furthermore we present the temperature dependent activation or deactivation of different Transient Receptor Potential (TRP) channels by means of planar patch clamping on our HTS platforms Patchliner and SyncroPatch 384PE as well as with highest resolution on a single channel level on our recently introduced Orbit mini setup.

CardioExcyte 96 Sensor Technology

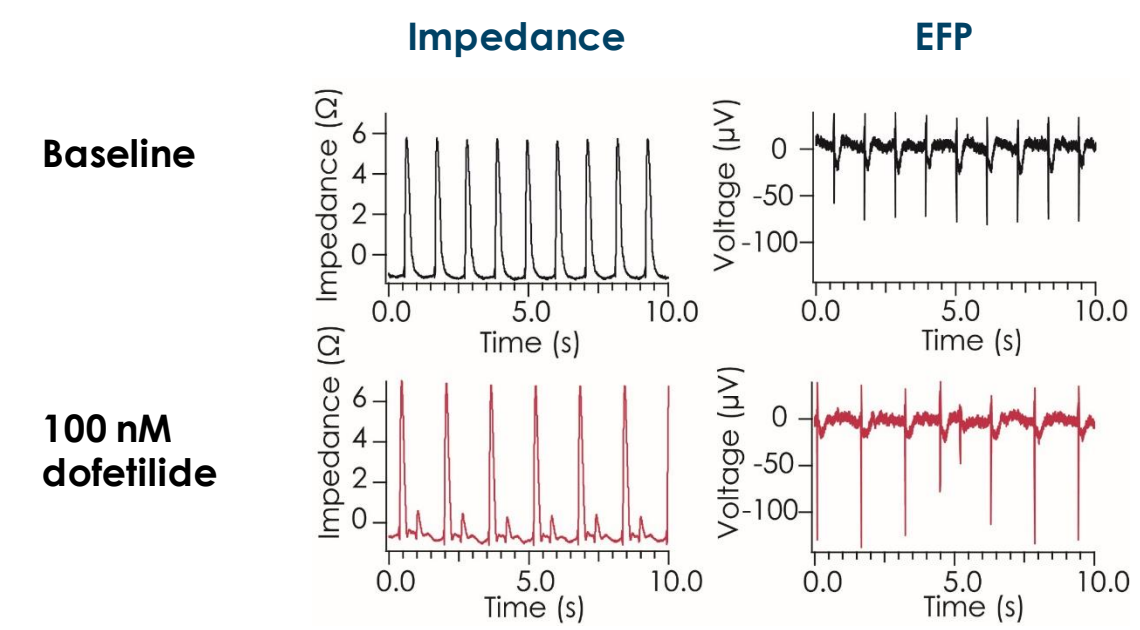


The CardioExcyte 96 combines impedance and EFP to measure effects on contractility and ion channel modulation (left). The NSP-96 plate contains gold electrodes embedded in each well (middle). The impedance (top) and EFP (bottom) signals are recorded from the same cell, an example is shown (right).

Development of the Impedance Signal



CardioExcyte 96 Compound Profiling



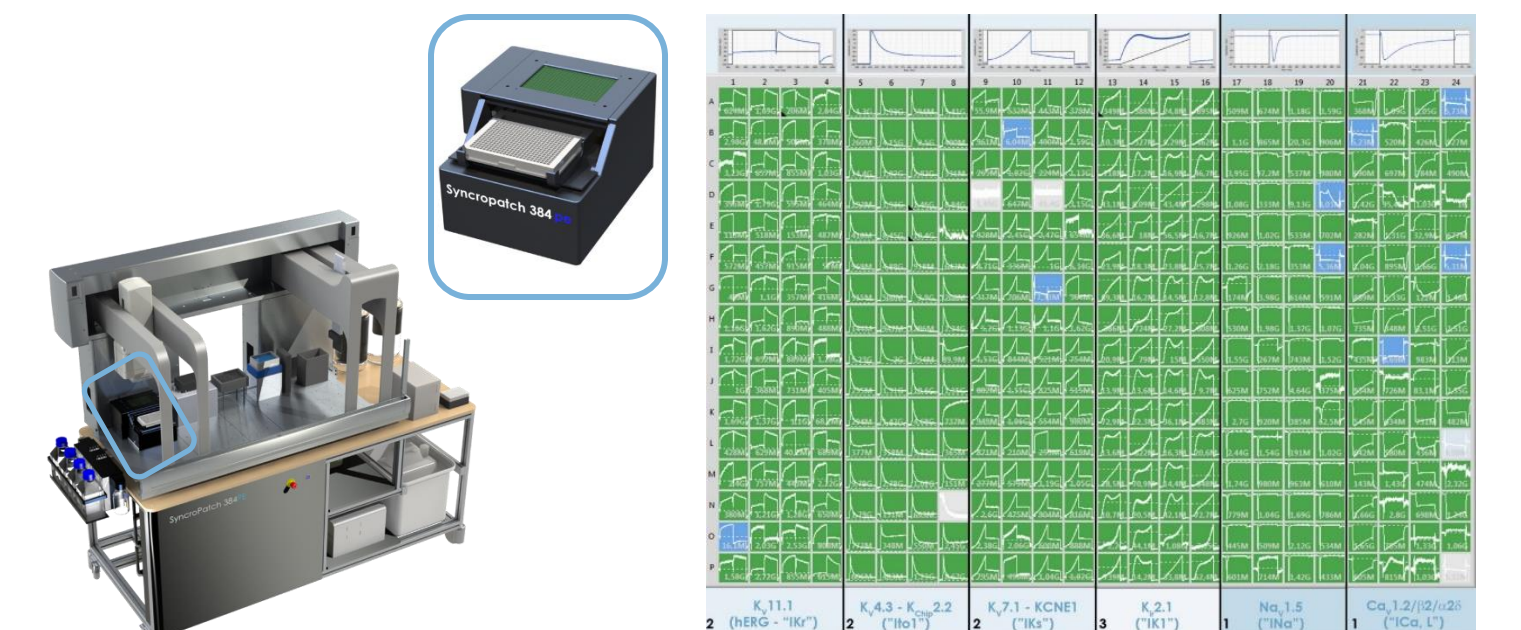
Dofetilide induces a decrease in beat rate and EADs in impedance (left) and EFP (right) modes.

Compound	Impedance			EFP			Effect
	AMP	Rate	RRD	AMP	Rate	RRD	
Dofetilide 100nM	+	-	-	+	-	-	ERG blocker
Cisapride 1 µM	+	-	-	+	-	-	ERG blocker
E-4031 1 µM	+	-	-	+	-	-	ERG blocker
Sotalol 1 µM	+	-	-	+	-	-	ERG blocker
Azelenolol 1 µM	+	-	-	+	-	-	ERG blocker
Guafendine 30 µM	+	-	-	+	-	-	ERG, K _v , Ca _v , Na _v blocker
Terfenadine 1 µM	+	-	-	+	-	-	ERG, Na _v blocker
Isoproterenol 30 µM	+	-	-	+	-	-	β-adrenergic agonist
Milrinone 300 µM	+	-	-	+	-	-	Ca _v blocker
Bay K 8644 1 nM	+	-	-	+	-	-	Ca _v activator

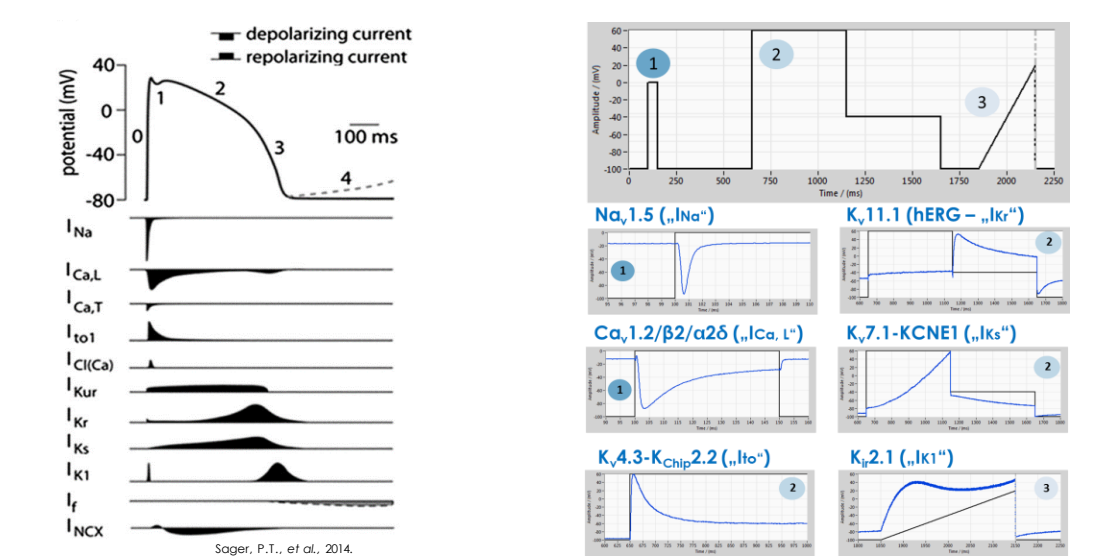
Ten different compounds were tested on Cor.4U cells in impedance and EFP modes. The results are summarized in this table.

Cor.4U Cardiomyocytes, AxioGenesis AG

Complementary Experiments Using the SyncroPatch 384PE



The SyncroPatch® 384PE is a patch clamp module that can be used with a state-of-the-art pipetting robot, e.g. Biomek FX (Beckman Coulter) shown here (left). Six different cardiac channels recorded simultaneously on the SyncroPatch 384PE. Screenshot shows data acquisition and analysis software performing recordings from HEK or CHO cells expressing the CiPA stipulated ion channels (ChanTest, a Charles River company) activated within one single experiment on the SyncroPatch 384PE.

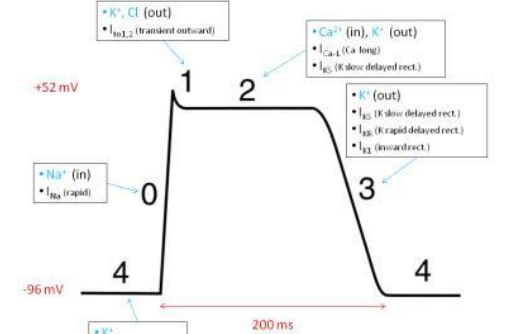


The ion channel currents mediating the cardiac myocyte action potential (AP; left). The voltage clamp protocol and examples of the 6 different currents recorded on the SyncroPatch 384PE.

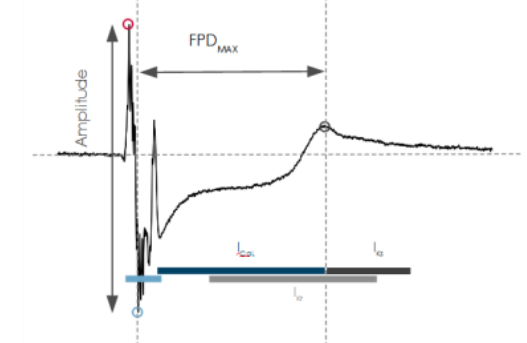
charles river

AP, EFP and ECG – How They Correlate

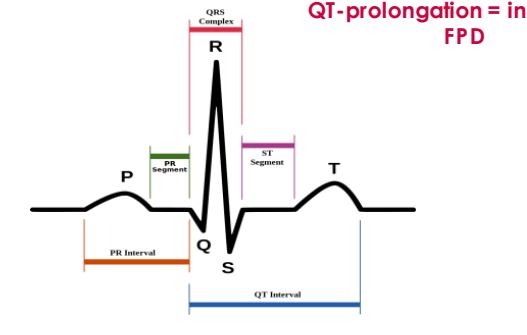
A. Human ventricular action potential¹



B. Extracellular electric field potential (EFP) and cardiac ion channel currents that contribute to iPSCM potentials.

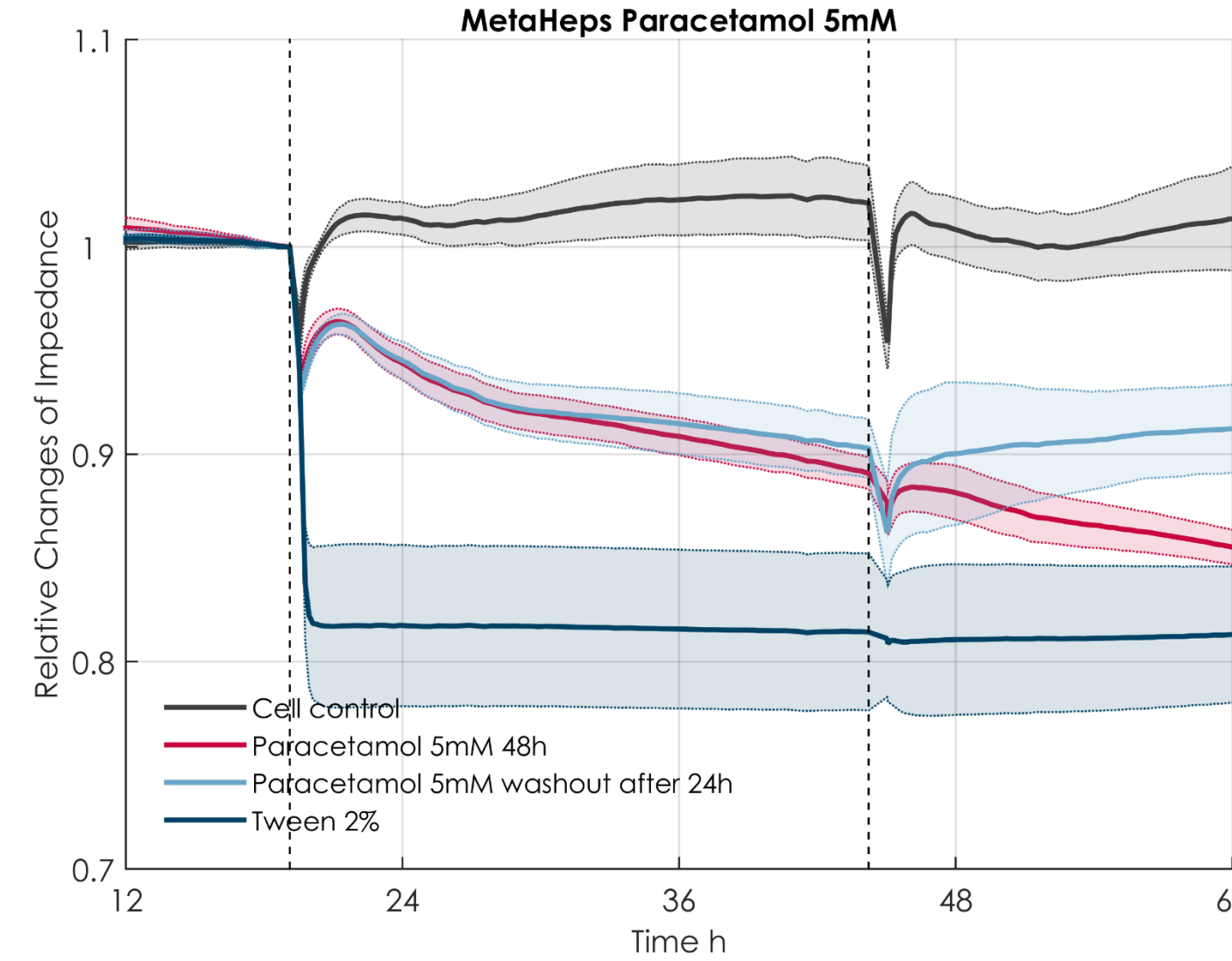


C. Surface electrocardiogram (ECG)



¹ Sherwood, L. (2022). Human Physiology: From Cells to Systems (9th (revised) ed.). Cengage Learning. ISBN 9781111577428.

Toxicology Screening on Hepatocyte-like Cells



Besides Cardiomyocytes, MetaHeps cells that exhibit hepatocyte-like characteristics have been studied with the CardioExcyte96. Application of 5 mM Paracetamol (APAP) led to a considerable decrease in overall impedance indicating severe cell damage within 24 hours of incubation (indicated by the dashed lines). Of note, removal of APAP after 24 hours resulted in detectable cell regeneration (light blue graph) while incubation with 5 mM APAP for 48 hours led to further decrease in impedance (pink graph).

MetaHeps®

Conclusions

- **CardioExcyte 96** is a hybrid system combining impedance and EFP recordings from the same cell.
- Recordings are made at physiological temperature in an incubator or using the Environmental Chamber.
- The **CardioExcyte 96** can be used in cardiac safety screening to measure effects of CiPA relevant compounds such as E4031.
- The **CardioExcyte 96** can be utilized for toxicology studies on e.g. hepatocytes or cell quality assessment.
- The **CardioExcyte 96** provides complementary data to other assays such as APC.



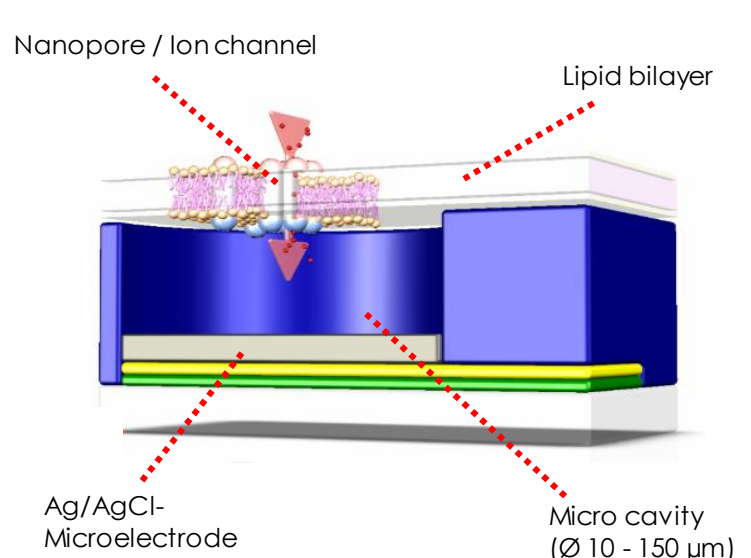
The Orbit mini with Temperature Control



- Including eFOUR 4 channel amplifier (Elements S.R.L., Cesena, I)
- Smallest footprint, USB powered
- Max. bandwidth: 100 kHz

elements
enabling technologies for Life Science

- Optional temperature control
- Active heating and cooling
- Temperature range : 5-50° C
- Software controlled



- Introducing unique MECA chip technology: micro electrode cavity array (lonera GmbH, Freiburg, D)
- Enabling recordings at highest bandwidth with excellent RMS noise behavior

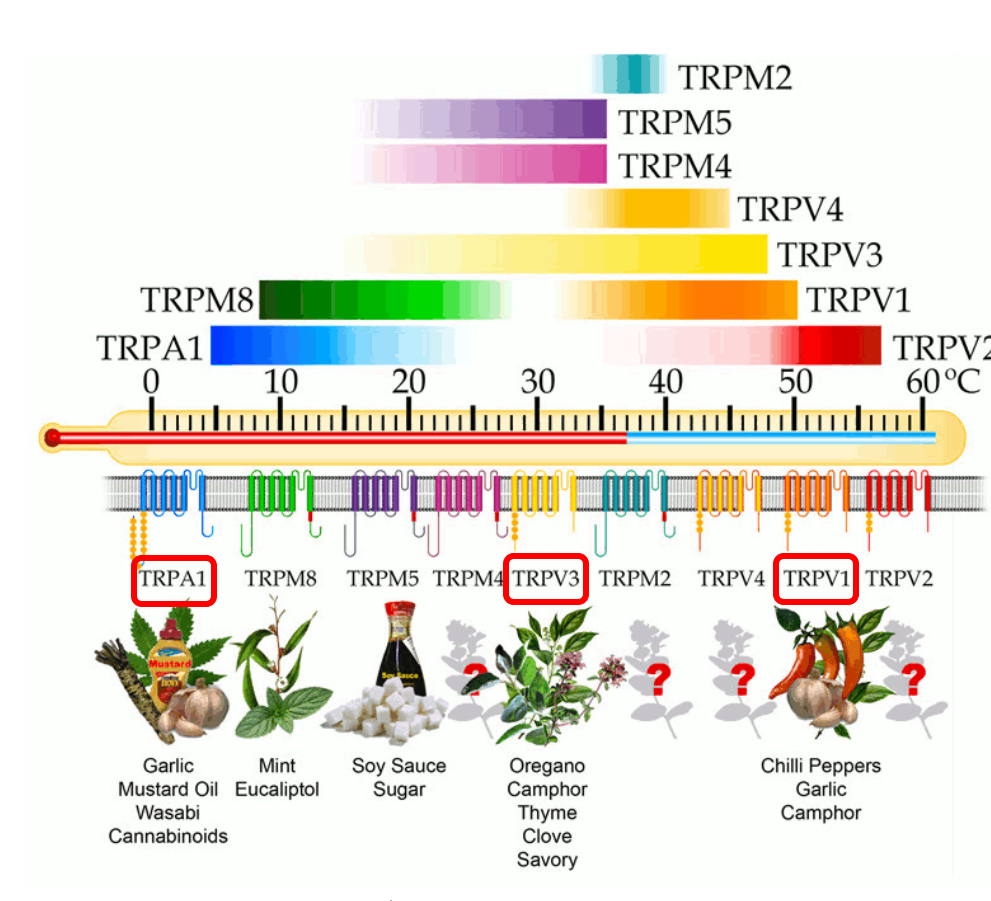
lonera

Typical Noise figures of a bilayer on:

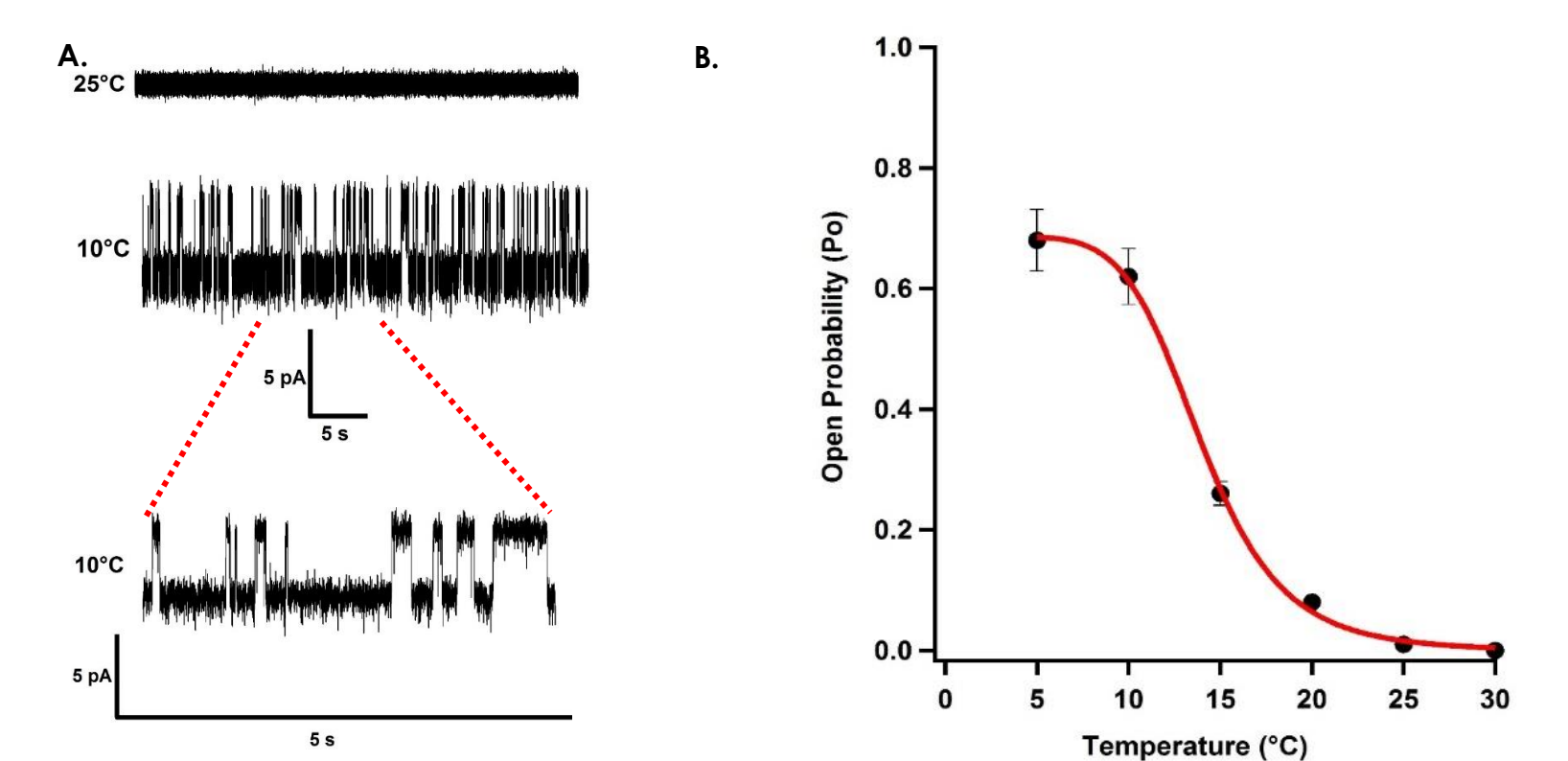
Orbit mini	Axon 200B amplifier
1.2 pA	950 fA – 5.5 pA (10 KHz)
7.0 pA	> 20 pA at (100 KHz) (depending on the bilayer setup used)

Thermo-sensitive TRP Channels

Thermal Transient Receptor Potential (TRP) channels are an important class of receptors found widely distributed throughout the mammalian central and peripheral nervous systems. They have been shown to be directly activated by heat or cold in physiologically relevant temperature ranges but are also activated by mechano-stimulation and various ligands. Understanding the mechanisms of temperature activation could lead to the discovery of novel compounds with differing effects on ligand activation and temperature activation for the treatment of pain and other disease states, with fewer side effects.

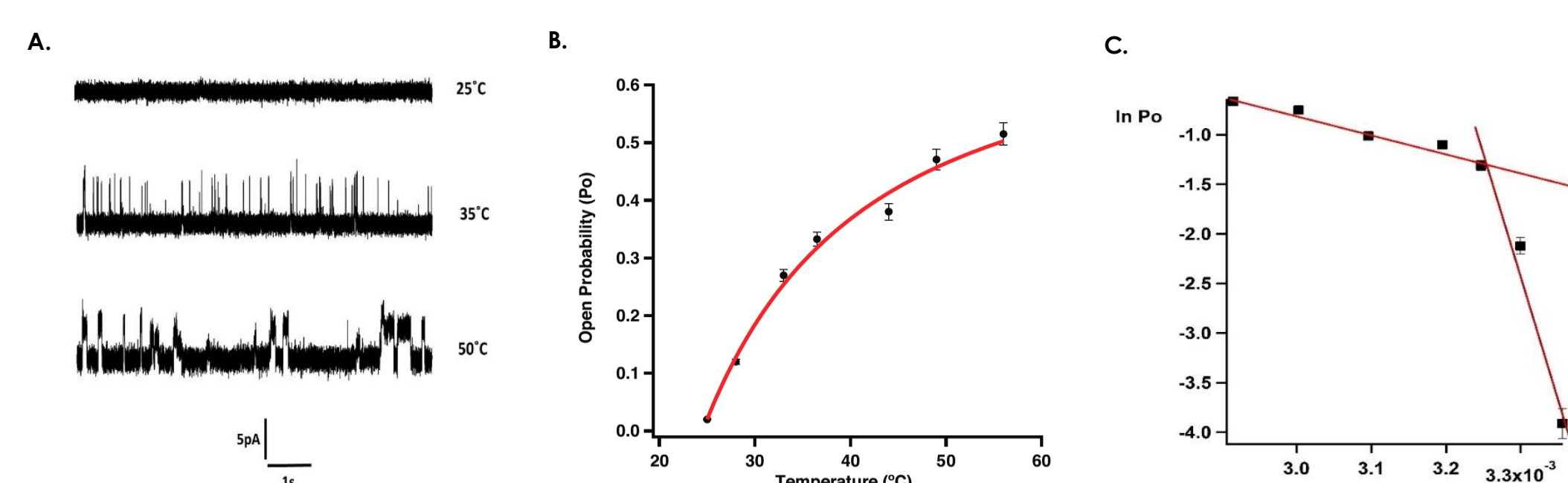


Cold Activated TRPA1



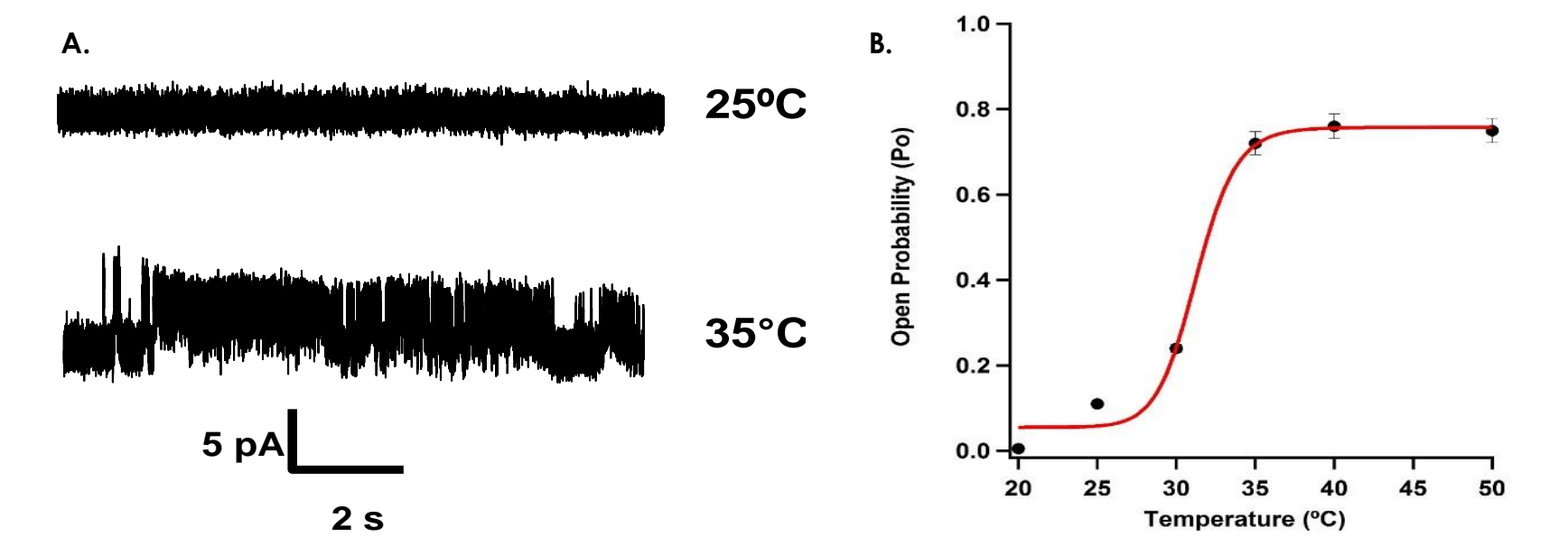
TRPA1 on the Orbit Mini. A. Effect of temperature on TRPA1 activity; B. The open probability (Po) versus the temperature and fitted with Boltzmann equation (EC₅₀ was found at 14°C). The Arrhenius plot of the same data resulted in a Q₁₀ of 46 (Literature: Q₁₀ ~ 40).

Heat Activated TRPV1



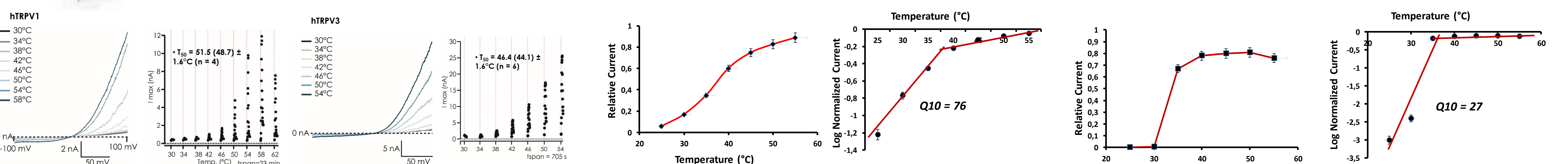
The effect of temperature on Purified TRPV1 reconstituted into lipid bilayers on the Orbit Mini. A. Evoking activity of TRPV1 in bilayers by elevated temperature; B. The effect of elevated temperatures on the open probability (Po) of TRPV1; C. Corresponding Arrhenius plot of the same data resulting in a Q₁₀ of 72 (Literature: Q₁₀ ~ 18 to >100).

Heat Activated TRPV3



Purified TRPV3 reconstituted into bilayers on the Orbit Mini. A. Effect of temperature on TRPV3 activity; B. Effect of elevated temperature on the open probability (Po). Fitting the data with the Boltzmann equation yields an EC₅₀ of 31.67°C. The Arrhenius plot of the same data results in a Q₁₀ of 22.2 (Literature: Q₁₀ ~ 17-56).

Correlation of Results Obtained via Automated Patch Clamp and via Bilayer Recordings with Temperature control



Response of TRPV1 and TRPV3 to an increasing temperature application in whole cell recordings on Nanion's Patchliner system. The T₅₀ values were determined at 51.5 °C and 46.4 °C respectively. Note that the overall conductance of the channels increases with the temperature.
Inducible TRPV1 CHO cell line kindly provided by Pfizer / TRPV3 HEK cell line kindly provided by Millipore

Thermodynamic analysis of TRPV1 relative current upon heat activation. Left: Current of TRPV1 obtained during a temperature increase. Right: temperature dependence presented in log(Po) versus temperature plot. The data resulted in a Q₁₀ of 76 (lipid bilayers experiments Q₁₀ = 72 and Literature: Q₁₀ ~ 18 to >100).

Thermodynamic analysis of TRPV3 relative current upon heat activation. Left: Current of TRPV3 obtained during a temperature increase. Right: temperature dependence presented in log(Po) versus temperature plot. The data resulted in a Q₁₀ of 27 (lipid bilayers experiments Q₁₀ = 22 and Literature: Q₁₀ ~ 17-56).

Literature values of TRP channels taken from 1. Mandrot, S., Roufogali, B.D. 2008. ThermoTRP Channels in Nociceptors: Taking a lead from Capsaicin receptor TRPV1. *Curr. Neuropharmacol.* 6:21-38. 2. Gavva, N.R., Treanor, J.J. et al. 2008. Pharmacological blockade of the vanilloid receptor TRPV1 elicits marked hyperthermia in humans. *Pain.* 136: 202-210. 3. Damien S.K. Sarwary and Terence M. Egan. Calcium dependent decrease in the single channel conductance of TRPV1. *Pflügers Arch.* 2011 Nov; 442(5): 681-691.