# Exploring stem cell-derived cardiomyocytes with automated patch clamp techniques

# RIKKE SCHRØDER<sup>1</sup>

# Abstract

There is increasing interest for cardiomyocytes as models for studying cardiac cellular physiology and preclinical drug safety testing. Stem cell-derived cardiomyocytes have the potential for such a model and have the possibility for modeling human diseases. The present investigation is the first to describe current properties from stem cell-derived cardiomyocytes using multi-hole recordings with planar automated patch clamp technology.

In our study pluripotent stem cell-derived cardiomyocytes were biophysically and pharmacologically characterized. The cells are differentiated in large numbers and cryo-preserved, which make them suitable for automated patch clamping and facilitate their use in drug screening. We tested the cells in two different recording modes; single-hole and multi-hole, respectively. For multi-hole recordings up to ten cells are patched at the same time and the total current is measured per site. This recording mode can be useful for small currents (e.g. endogenous) and typically increases the success rate for useful data. For all experiments the whole-cell configuration was used and three different types o currents were studied; Na<sup>+</sup>, Ca<sup>2+</sup> and K<sup>+</sup>. Using specific voltage protocols biophysical characteristics of each current was described and compared from single-hole and multi-hole experiments. We showed that currents recorded from these pluripotent stem cell-derived cardiomyocytes are similar to human cardiomyocytes and the response to known pharmacology is as expected. The  $V_{0.5}$  values, I-V related onships, current kinetics and IC<sub>ro</sub> values determined for known blockers (TTX, nifedipine and cisaprid</sub>were comparable for the two recording modes. Clearly the success rate for usable data per measu rement plate was significantly increased with the multi-hole technology. This is the first time current properties of stem cell-derived cardiomyocytes have been described from multi-hole recordings with planar automated patch clamp. Our study has shown that automated patch clamp is ready for stem cell-derived exploration.

### Material and Methods

#### **Ringer solutions voltage clamp experiments** EC I<sub>Na</sub> (mM): 120 NaCl, 5 KCl, 3.6 CaCl<sub>2</sub>, 1 MgCl<sub>2</sub>, 20 TEACl, 10 HEPES. pH 7.4 with NaOH nM): 15 NaCl, 140 KCl, 1.2 CaCl<sub>2</sub>, 1 MgCl<sub>2</sub>, 10 HEPES. pH 7.4 with NaOH EC I<sub>ca</sub> (mM): 125 NMG, 3.6 CaCl<sub>2</sub>, 1 MgCl<sub>2</sub>, 20 TEA Chloride, 10 HEPES. pH 7.4 with HCl

IC I<sub>ca</sub> and I<sub>Na</sub> (mM): 120 CsCl, 3 MgCl<sub>2</sub>, 10 EGTA, 5 HEPES, 4 Na<sub>2</sub>-ATP. pH 7.4 with CsOH IC for for  $I_{\kappa}$  (mM): 5.374 CaCl<sub>2</sub>, 1.75 MgCl<sub>2</sub>, 3.125/10 KOH/EGTÁ, 120 KCl. pH 7.2 with KOH

Ringer solutions current clamp experiments EC (mM): 145 NaCl, 4 KCl, 2 CaCl, 1 MgCl, 20 TEACl, 10 HEPES, 10 glucose. pH 7.4 with NaOH IC (mM): 5.374 CaCl<sub>2</sub>, 1.75 MgCl<sub>2</sub>, 3.125/10 KOH/EGTA, 10 HEPES, 120 KCl, 4 Na<sub>2</sub>-ATP. pH 7.2 with KOH

# Cells

Cryopreserved human induced pluripotent stem (iPS) cell-derived cardiomyocytes, iCells<sup>®</sup>, were generously received from Cellular Dynamics. Each vial contained approximately 4.5 mio cells with a plating efficiency of 41-53%. The cells were thawed and plated on fibronectin coated T-25 culture flasks using 60,000 cells per cm<sup>2</sup> to obtain a syncytial monolayer. The culture flasks were stored in a controlled environment in an incubator (37°C, and 5%  $CO_2$ ) and the culture media was exchanged every other day. The culturing was continued for six to eight days. After culturing the cells were harvested and prepared for QPatch experiments according to the SOP developed by Sophion.

For action potential recordings human embryonic stem (hES) cell-derived Cytiva™ Cardiomyocytes were generously received from GE Healthcare. Cells were received in frozen vials. Each vial contained approximately 2 mio cells with a purity of 64%. The cells were thawed and plated on fibronectin coate T-25 culture flasks using 60.000 cells per cm<sup>2</sup> to obtain a syncytial monolayer. Cells were cultured according to the SOP received from GE Healthcare. After culturing the cells were harvested and prepared for QPatch experiments according to the SOP developed by Sophion.

### QPatch

Small volumes of 50-100 µl cell suspension were applied to the QPatch. Cell positioning, giga sealing and whole cell formation were obtained automatically with suction pulses according to a pre-program med suction protocol. Currents were recorded using either the single-hole or multi-hole technology. For the multi-hole technology up to 10 cells are patch per measurement site and the total current is recorded. For the single-hole technology one cell is patched per measurement site. Voltage protocols were applied to obtain specific biophysical characteristics of the currents. One or more concentrations of the compounds were added sequentially and automatically through the flow channels in the QPlate. Action potential recordings were obtained after giga seal and whole cell formation. Data were analyzed using the QPatch Assay Software.





### **References:**

Schneider M, Proebstle T, Hombach A, Rüdel R Characterization of the sodium currents in isolated human cardiocytes. Plügers Arch, 1994;84-90

Drolet B, Khalifa M, Daleau P, Hamelin B, Turgeon J Block of the rapid component of the delayed rectifier potassium current by the prokinetic agent cisapride underlies drug-related lengthening of the QT interval. Circulation, 2011;204-210

Sanguinetti MC, Jurkiewicz NK Two components of cardiac delayed rectifier K+ current. J Gen Physiol, 1990;195-215

Pelzmann B, Schaffer P, Bernhart E, Lang P, Mächler H, Rigler B, Kiodl B L-type calcium current in human ventricular myocytes at a physiological temperature from children with tetralogy of fallot. Cardiovascular Research, 1998;424-432

GE Healthcare Cell Technologies The Maynard Centre, Whitchurch Cardiff CF14 7YT UK

MORTEN SUNESEN<sup>1</sup>





I-T plot;  $I_{Nav}$  current blocked by increasing concentrations of TTX (1, 3, 10, 30  $\mu$ M)





Boltzmann fit of  $I_{Na}$  tail current obtained at +10 mV from currents pre-stimulated with voltage steps from -100 mV to +70 mV. The  $V_{0.5}$  was -49.4 mV (± 2.9, n=5). (V<sub>0.5</sub> lit. value -62 to -72 mV, Schneider et al. 1994)<sup>33</sup>



I-T plot; I<sub>c</sub>, current blocked by increasing concentrations of nifedipine (0.1, 1, 10



4000 0 60 100 160 200 260 300 350 -50 -40 -30 -20 -10 0 10 20 30 40 50 Step Voltage [mV] 746.1.1'4:[0]Break-in ♥ Min(Ca-act) I-V plot of the  $I_{c_2}$  current.  $I_{c_2}$  was activated by voltage steps from -60 mV to +50 mV. The current began to activate close to -40 mV and peaked at voltages between

0 and +10 mV. (Similar I-V relations, Pelzmann et al. 1998)



The study of ion channels in their native tissue using automated patch clamp has historically been impeded by factors like; the cells have not be available in sufficient numbers and the cell population has shown inadequate homogeneity. With the recent efforts in deriving cardiomyocytes from pluripotent-, and embryonic stem cells, cardiomyocytes are now available in adequate quantities and with acceptable purity to match a range of applications for automated patch clamp technology. Still, the stem cell-derived cardiomyocytes exhibit some variation in expression of specific current sub-types, which limits the data point throughput for automated patch clamp systems.

Here we show single-hole and multi-hole recordings of I<sub>Na</sub>, I<sub>Ca</sub>, I<sub>K</sub> recorded from human iPS cardiomyocytes. For the first time iCell cardiomyocytes have been recorded using the multi-hole technology for automated patch clamping. Our results showed that biophysical and pharmacological characteristics of I<sub>Na</sub> were comparable using the single-hole and the multi-hole technique, respectively. The success rate for experiments with useable I<sub>Na</sub> significantly increased using the multi-hole technology. From our single-hole recordings we found that most cells expressed I<sub>Na</sub> and  $I_{\kappa}$ , and less than half of the cells expressed  $I_{Ca}$ . The recordings from  $I_{Na}$   $I_{Ca}$ ,  $I_{\kappa}$  were similar to recordings obtained and reported from other mammalian cardiomyocytes.

We used hES Cytiva cells for current clamp recordings obtained on the QPatch system. These cells exhibited spontaneous action potentials with ectrophysiological properties expected for ventricular subtypes (similar data obtained from iPS iCell cardiomyocytes, data not shown).

In conclusion, stem cell-derived cardiomyocytes are commercially available in large quantities, are easy to handle and show satisfactory success rates especially for multi-hole recordings with the QPatch automated patch clamp system. The iPS and hES cells used in this study have shown expected biophysical and pharmacological behavior from voltage clamp and current clamp recordings. From our exploration of different stem cell-derived cardiomyocytes our data have shown that these cells are candidates for in vitro cardiac electrophysiological studies and can be a power full tool for cardiac safety research.



**CELLUIAR Dynamics** international

# CELLULAR DYNAMICS<sup>2</sup> INTERNATIONAL, INC. 525 Science Drive Madison, WI 53711 UNITED STATES





iCell cardiomyocytes cultured for 7 days in a fibronectin coated culture flask



QPlate sealing overview from a **single-hole** experiment. The success rate

for sealing for single-hole experiments could reach up to 70 % for sealing

and 60 % for whole-cells

iCell cardiomyocytes harvested by trypsination procedure. The cell density is approximately 1.5 mio cells per ml.

	Single-hole	
Seal	53 % (±12, n=9)	
Whole-cell	22 % (±13, n=9)	

Table 1. Success rates for obtained seals and whole-cells with iCells n refers to the number of QPlates 16's used for single-hole experiments

	Single-hole	Multi-hole
Cell size	29 pF (±15, n=41)	N.D
Cells expressing I <sub>Na</sub>	91 % (±16, n=40)*	N.D
Cells expressing I <sub>ca</sub>	41 % (±39, n=58)*	N.D
Cells expressing I <sub>k</sub>	80 % (±35, n=17)*	N.D
Usable I <sub>Na</sub> data/QPlate	16 % (±10, n=224)	58 % (±22, n=80)
I <sub>Nav</sub> amplitude	-3.5 nA (±2.2, n=24)	-6.4 nA (±42., n=46)
IC <sub>50</sub> TTX for I <sub>Na</sub>	10.3 µM (±0.5, n=6)	6.3 µM (±4.5, n=12)
Tau for I <sub>Na</sub> inactivation***	0.85 ms (±0.29, n=22)	0.99 ms (±0.16, n=20)

#### Table 2. Comparing different re sults from $I_{Na}$ , $I_{Ca}$ and $I_{\kappa}$ obtained from single-hole and multi-hole recordings of iCell Cardiomyocyte \*Specific currents from established whole-cells. \*\*Recorded at 0 mV. \*\*\*Data fitted to a monoexponentia curve



Spontaneous action potential recorded from Cytiva<sup>™</sup> Cardiomyocytes. A beating frequency of approximately 1 Hz was observed. Each action potential had an amplitude of 80 mV and duration of 400 ms

# Conclusion