



Application Report

hiPSC-derived cardiomyocytes: A robust and reproducible QPatch® II assay for Na_v1.5 measurements

Including biophysical characterization and concentration-dependent block of Na_v1.5

Summary

The Na_v1.5 channel is critical in maintaining the physiological electrical function in cardiomyocytes (CMs). Consequently, high-throughput measurements of Na_v1.5 currents (I_{Na}) in hiPSC-CMs are useful e.g. in the characterization of cardiac disease phenotypes and compound testing. QPatch II has been demonstrated as a capable platform for such measurements¹, and here we present a high-performance Na_v1.5 assay using iCell® Cardiomyocytes² (hiPSC-CMs) from FUJIFILM Cellular Dynamics with:

- Assay success rates ~45% with more than 85% hiPSC CMs expressing I_{Na} with peak currents $|I_{Na}| > 200$ pA (at -30 mV)
- Biophysical characterization with activation/inactivation potentials in agreement with literature values
- Cumulative compound concentration-response of the local anesthetic agent tetracaine with IC₅₀ values in agreement with literature values

Results and discussion

Assay success rate and measurement quality

We obtain an assay success rate of ~45%, calculated as the percentage of measurement sites, out of the full measurement plate (48 sites), that pass our quality criteria (membrane resistance, $R_{mem} > 200$ M Ω , cell capacitance, $C_{slow} > 4$ pF). In 85% of hiPSC-CMs we record I_{Na} with peak current $|I_{Na}| > 200$ pA. See Figure 1 for a partial measurement plate view, displaying I_{Na} current-voltage (IV) relationship recorded in 42% of the sites.

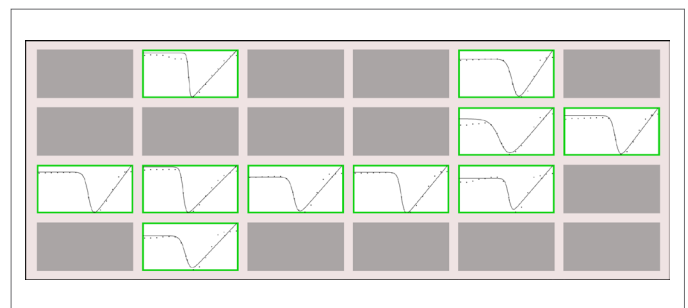


Fig. 1: Plate overview of half a QPatch consumable (24 measurement sites out of 48) displaying I_{Na} IV relationship recorded in individual hiPSC-CMs. In this experiment we obtained a 42% success rate.

R_{mem} and C_{slow} values are plotted in Figure 2A for all recorded hiPSC-CMs (~30% of 48 sites) that maintained a good seal for the duration of the concentration-response experiment. Note that hiPSC-CMs display a large distribution of sizes as observed in the C_{slow} values that varied from 4 pF to 40 pF (Fig. 2A, right). To ensure good voltage clamping we reduced the extracellular concentration of Na⁺, see Figure 2B for a representative I_{Na} trace. The peak I_{Na} was recorded at -30 mV and correlated well with cell size. See Figure 2C for plots of Na_v1.5 peak current, I_{Na} , and peak current density.

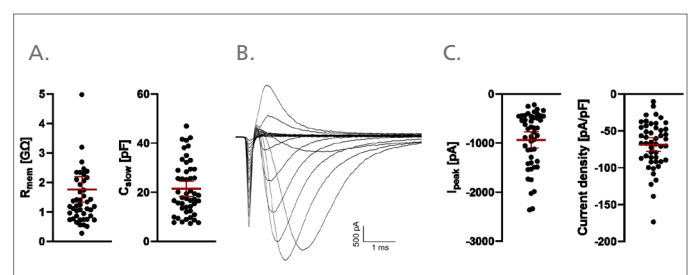


Fig. 2: Measurement quality and representative hiPSC-CM I_{Na} current. **A)** Seal resistance, R_{mem} [G Ω], (left) and cell capacitance, C_{slow} [pF], (right) (mean and 95% CI are shown in red, $N_{CMs} = 40$, $N_{QPlates} = 3$). **B)** Representative Na_v1.5 current trace in response to a voltage step protocol (see methods). **C)** Na_v1.5 peak current, I_{Na} [pA], (left) and peak current density [pA/pF] (right) (mean and 95% CI are shown in red, $N_{CMs} = 48$, $N_{QPlates} = 3$).

Biophysical characterization

We quantified the biophysical properties of the I_{Na} recorded from hiPSC-CMs. The experiments were conducted at a reduced extracellular sodium concentration ($[Na^+] = 10$ mM), to ensure good voltage clamping of the cells. The I_{Na} activated around -60 mV and peaked around -30 mV (Fig. 3A). The half maximal activation was ~ -44 mV and the half maximal inactivation was ~ -83 mV, in agreement with reported values for $Na_v1.5^{2,3}$ (Fig. 3B).

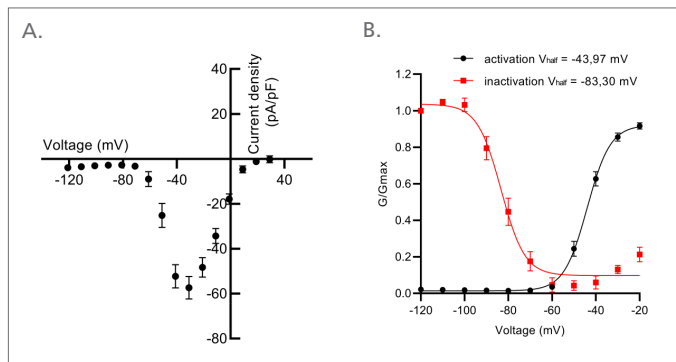


Fig. 3: Biophysical characterization of the I_{Na} current. **A)** Average I_{Na} current density versus step voltage (voltage steps of 10 mV from a holding potential of -130 mV, see Methods) ($N_{CMs} = 40$). **B)** Normalized conductance (G/G_{max}) versus step voltage was plotted for the activation (black, $N_{CMs} = 39$) and inactivation (red, $N_{CMs} = 49$) and fitted with the Boltzmann equation (solid lines) to yield the voltage at half maximum activation ($V_{half} \sim -44$ mV) and inactivation ($V_{half} \sim -83$ mV). All data points are mean \pm sem.

Cumulative concentration-response of tetracaine

As a proof-of-concept concentration-response experiment, we recorded the I_{Na} sensitivity to increasing concentrations (0.01 μ M to 100 μ M) of the local anesthetic agent tetracaine, using a two-pulse voltage protocol, including an adaptive protocol for V_{half} measurement for improved control of $Na_v1.5$ state before measurement⁴ (see Methods). Tetracaine blocked I_{Na} in a state-dependent manner, and in agreement with previous APC measurements on immortal cell lines expressing the $Na_v1.5$ ion channel⁵, we quantify an IC_{50} value of ~ 2.4 μ M for the second peak of the double pulse.

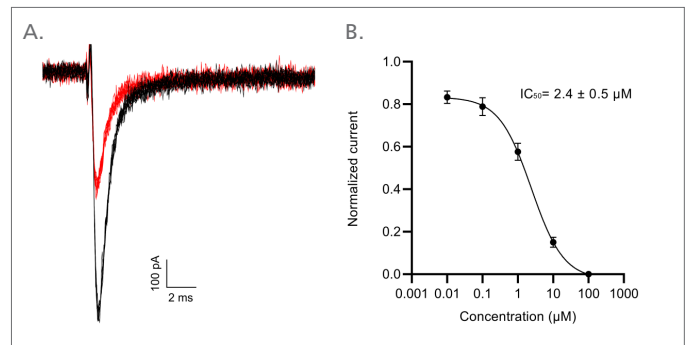


Fig. 4: Concentration-dependent block of I_{Na} in hiPSC-CMs. **A)** Representative I_{Na} current (recorded in the second pulse of the voltage protocol) before (black) and after (red) the addition of 1 μ M tetracaine. **B)** Normalized current (I_{Na}) as a function of tetracaine concentration ($N_{CMs} = 24$). Fitting of the Hill equation (solid line) to the data yielded an IC_{50} value of ~ 2.4 μ M. Data points are mean \pm sem.

Results and discussion

In conclusion, we demonstrate a successful $Na_v1.5$ assay using iCell Cardiomyoctes² (hiPSC-CM) from FUJIFILM Cellular Dynamics, with assay success rates up to 45%. The cell line shows robust and reproducible I_{Na} current in more than 85% of recorded hiPSC-CMs. The biophysical and pharmacological features measured using QPatch II, are in concordance with I_{Na} previously reported, from hiPSC-CMs^{3,6}.

Methods

iCell Cardiomyocytes² (hiPSC-CMs) were kindly provided by FUJIFILM Cellular Dynamics. They were cultured according to vendor SOP and assayed (7 – 21) days post-thawing the frozen vials. The Na_v1.5 current density was stable within this time frame.

Cell dissociation

On experiment day the cells were dissociated using a Sophion SOP, specifically developed for iPSC-CMs and further optimized for iCell Cardiomyocytes². For more information contact the authors directly or through Sophion at info@sophion.com.

Solutions

Extracellular solution (in mM): 2 CaCl₂, 1 MgCl₂, 10 HEPES, 4 KCl, 10 NaCl, 135 C₅H₁₄NINO, 10 Glucose, pH = 7.4 with NaOH, osmolarity = 305 mOsm with sucrose

Intracellular solution (in mM): 10 NaCl, 10 HEPES, 1 EGTA, 5 CsOH, 140 CsF, pH = 7.3 with HCl, osmolarity = 280-290 mOsm with sucrose

Whole-cell protocol

We used a whole-cell protocol, specifically developed for iPSC-CMs. For more information contact the authors directly or through Sophion at info@sophion.com.

Voltage protocols

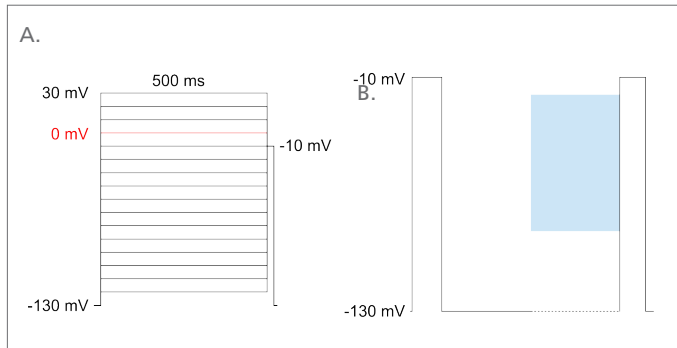


Fig. 5: Applied voltage protocols. A) Voltage step protocol for recording I_{Na} current-voltage relationship. The conductance (G) was calculated as $G = I / (V - V_{rev})$, with the activation current being measured during the 500 ms step pulses from -120 mV to +30 mV, and the steady state fast inactivation current being measured during a 20 ms test pulse to -10 mV following the pre-pulse step voltages. B) Double pulse for measurement of state-dependent I_{Na} compound effects. Individually recorded V_{half} values (voltage at half maximum activity) were applied before the second pulse (illustrated by the blue square) using our adaptive protocol for measurements of V_{half}⁴.

References

1. Badone, B. & Karatsiompani, S. hiPSC-derived cardiomyocytes - Nav1.5 compound screening on QPatch® II. <https://shorturl.at/agolM>. (2022).
2. Moreau, A. et al. Sodium overload due to a persistent current that attenuates the arrhythmogenic potential of a novel LQT3 mutation. *Front Pharmacol* 4 OCT, (2013).
3. Lemoine, M. D. et al. Human Induced Pluripotent Stem Cell-Derived Engineered Heart Tissue as a Sensitive Test System for QT Prolongation and Arrhythmic Triggers. *Circ Arrhythm Electrophysiol* 11, (2018).
4. Kammonen, J. & Korsgaard, M. P. G. VHalf adaptive protocols on Qube. <https://shorturl.at/btJWZ> (2019).
5. Sauter, D., Christensen, M. T., Jacobsen, R. B. & Mattsson, G. Biophysical and pharmacological profiling of multiple voltage-gated sodium channel subtypes on QPatch II. Poster presented at Ion Channel Modulation Symposium Preprint at <https://shorturl.at/wMRW4> (2019).
6. Van De Sande, D. V. et al. Pharmacological profile of the sodium current in human stem cell-derived cardiomyocytes compares to heterologous Nav1.5+β1 model. *Front Pharmacol* 10, (2019).



iCell Cardiomyocytes² (CM²) were kindly provided by FUJIFILM Cellular Dynamics, Inc.

www.fujifilmcdi.com

Authors:

Kadla Røskva Rosholm, Senior Research Scientist
Stefania Karatsiompani, Application Scientist
Ozhathil Lijo Cherian, Research Scientist R&D