

Application Report

HEK-hKir2.1 tested on QPatch

The biophysical characteristics of the hKir2.1 channel were studied from high resistance whole cell recordings in IV- and dose-response experiments and the values correspond to literature values.

Introduction

The aim of this report is to demonstrate the performance of HEK-hKir2.1 from CreaCell on QPatch. The cell line was characterized in terms of biophysical properties such as IV-relationship, effect of extracellular K^+ , effect of Ba^{2+} and pharmacologically properties of the channel.

Results and discussion

Experiments were carried out using four different protocols in order to study 1) IV-relationship, 2) effect of extracellular K^+ and 3) effect of the modulator chloroethylclonidine (CEC).

The IV-relationships for was studied by applying a step protocol from -150 mV to 0 mV with a holding potential at -20 mV, where the hKir2.1 channels are non-conducting. Stepping to potential negative to the reversal potential result in a large inward current. At potential more positive than E_{rev} result in a small outward current. Raw current traces from the current-voltage relationship experiment are shown in figure 1a. The corresponding IV curve is shown in figure 1b. The reversal potential is observed at $E_{rev} = -80.2 \pm 0.54$ mV, $n=8$ which corresponds to the theoretical E_{rev} that is -85.9 mV for this channel under these specific experimental conditions. The maximal inward current response $I_{max} = -2824.23 \pm 234.24$ pA, $n=8$ was observed at a potential of $V_{min} = -138.15 \pm 1.96$ mV, $n=8$. The small inward current response was observed at $I_{min} = 143.81 \pm 12.01$ pA, $n=8$ at a potential of $V_{max} = -62.16 \pm 0.99$ mV, $n=8$.

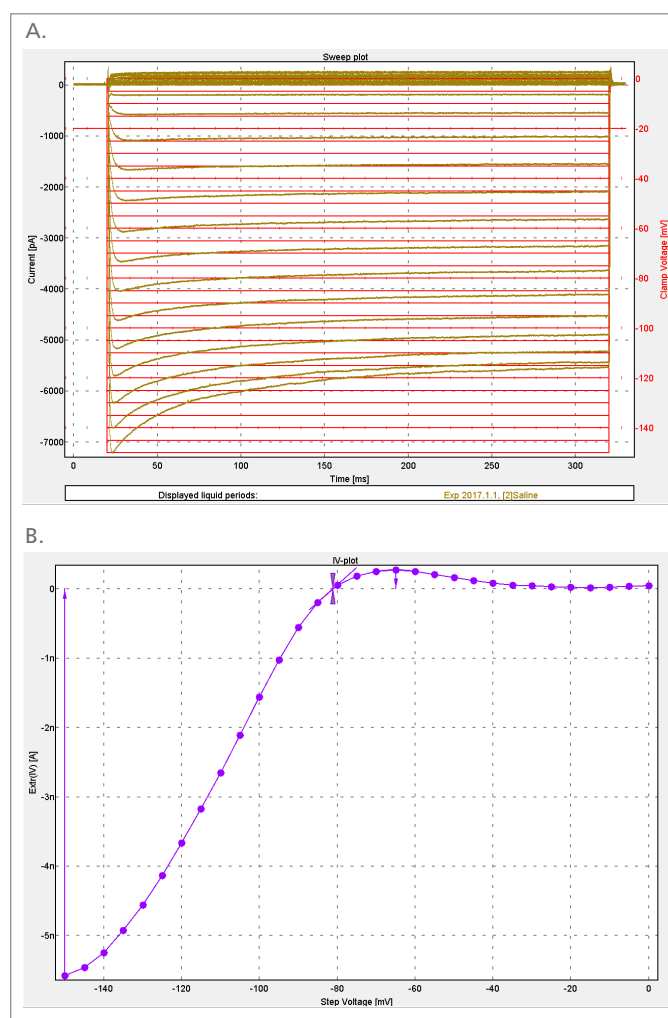


Fig. 1 : A. Raw hKir2.1 current traces during a step voltage protocol from -150 mV to 0 mV with 5 mV step increment. B. IV curve showing a strong inward current a low potential and a small outward current at potentials positive to E_{rev} .

Effect of external K⁺

In the next section, we show the effect of increasing K⁺ in the extracellular solution. By increasing the K⁺-concentration from 4 mM to 8 and 16 mM K⁺ in the extracellular solution, an increase in the conductance is observed in both the inward (negative to E_{rev}) and the outward (positive to E_{rev}) direction in hKir2.1 (Figure 2). At the same time, a parallel change in E_{rev} was observed in accordance to N_{ernst} equation for 4 mM K⁺=-84.2 mV, 8 mM K⁺=-69.0 and 16 mM K⁺=-55.7 mV (theoretical; 4 mM K⁺=-85.9 mV, 8 mM K⁺=-68.4 mV and 16 mM K⁺=-50.9 mV, 1).

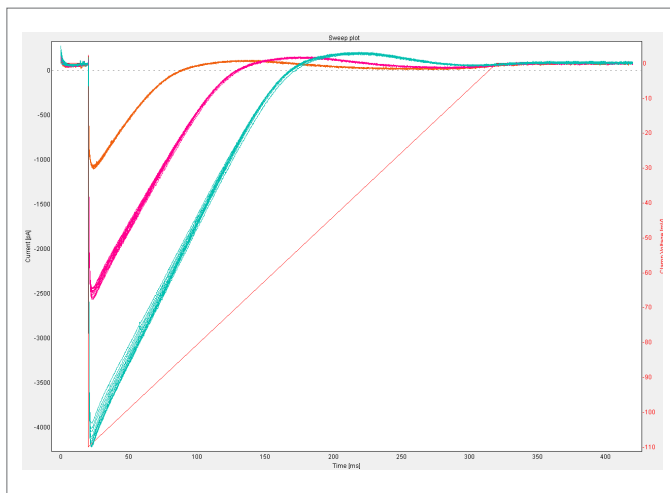


Fig. 2: Effect of extracellular K⁺. Raw current traces at 4 (orange), 8 (pink) and 16 mM (blue) external K⁺.

Effect of Ba²⁺ on hKir2.1

The hKir2.1 channel is blocked by external Ba²⁺ in a voltage-dependent manner. Figure 3a shows the effect of 5 increasing concentrations of external Ba²⁺, starting from 100 mM in a 10-fold dilution. The corresponding Hill fit is shown in Figure 3b. IC₅₀ is estimated to 5.12±0.42 μM, n=10, which corresponds well to literature values (2).

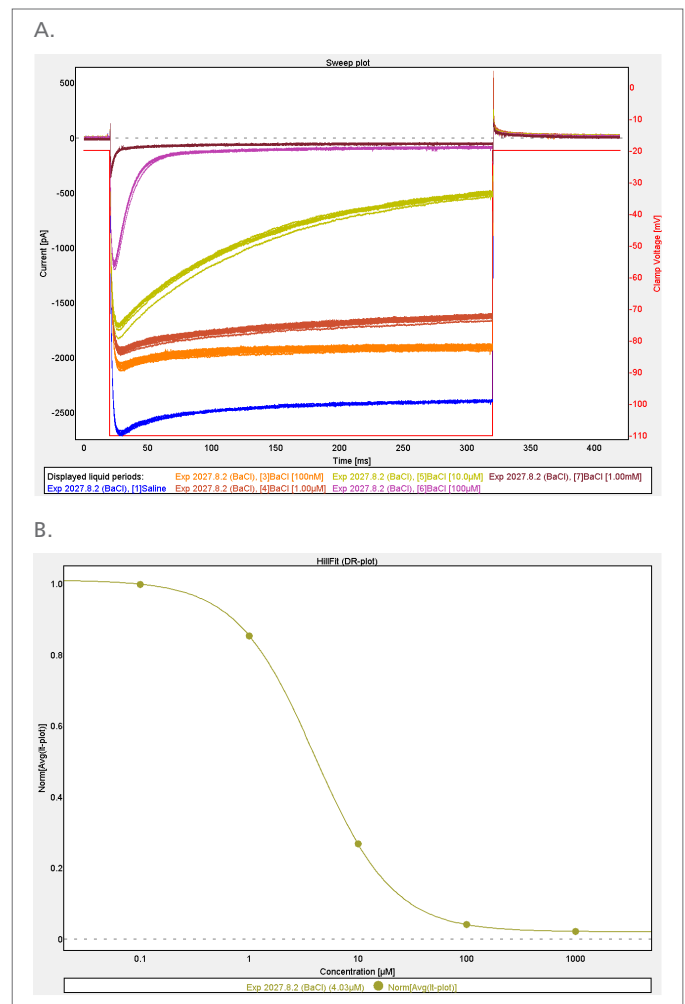


Fig. 3: A. Raw current traces from application of 5 increasing concentrations of extracellular Ba²⁺. B. Normalized Hill fit showing the dose-dependency of Ba²⁺ on hKir2.1.

Pharmacological effect of Chloroethylclonidine (CEC)

In the next section we will show data from pharmacological voltage-dependent block of hKir2.1 by CEC. The experiment was performed by application of extracellular solution with four increasing concentrations of CEC: 15.6 μM, 63 μM, 250 μM and 1 mM. Raw current traces are shown in Figure 4.

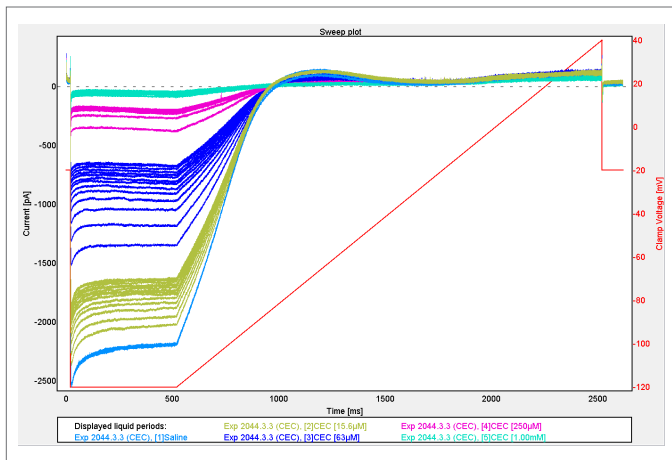


Fig. 4: Raw current traces of hKir2.1 with four increasing concentrations of CEC applied.

Figure 5a shows the time-current plot and Figure 5b the corresponding Hill fit. In average the IC_{50} for CEC was estimated to $IC_{50}=31.03\pm 0.97$, $n=11$.

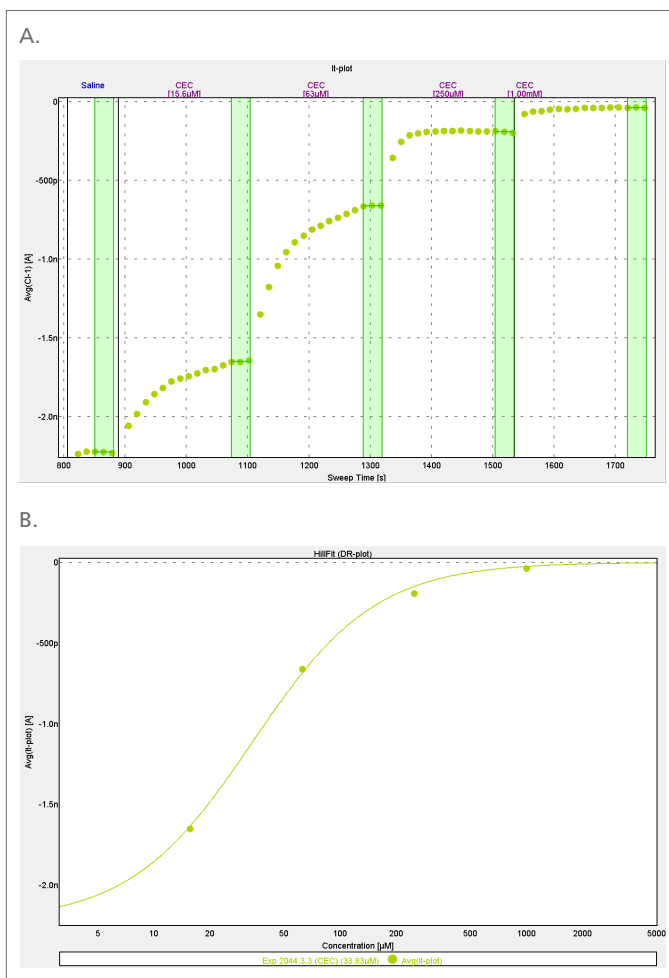


Fig. 5: A. Time-current plot from a single cell with four increasing concentrations of CEC.. B. Hill fit showing four concentrations of CEC.

Conclusion

We have demonstrated the functionality of HEK-hKir2.1 on the QPatch. Biophysical characteristics of the hKir2.1 channels were studied from high resistance whole cell recordings in IV- and dose-response experiments and the values found on the QPatch correspond well to published literature values.

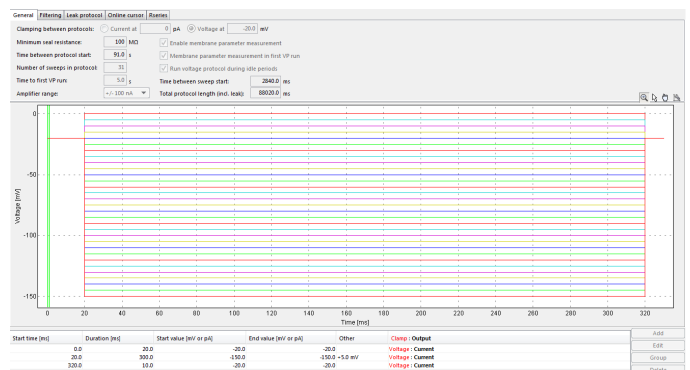
Methods

HEK-hKir2.1 cells were grown and harvested according to the SOP's specified from CreaCell (CreaCell.com) and modified for use on QPatch by Sophion Bioscience.

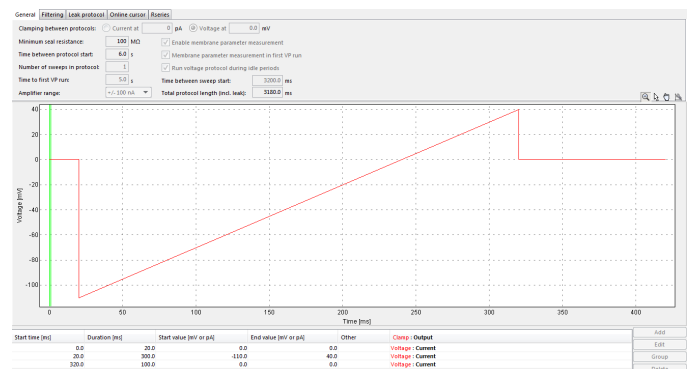
Voltage protocols

For experiments with HEK-hKir2.1 the following protocols were used. 1) Step IV protocol from -150 mV to 0 mV in 5 mV steps, $V_{hold} = -20$ mV, 2) Ramp IV protocol from -110 mV to +40 mV, $V_{hold} = 0$ mV, 3) short pulse from -20 mV to -110 mV for 300 ms, $V_{hold} = -20$ mV and 4) Long ramp protocol with a 500 ms step at -120 mV.

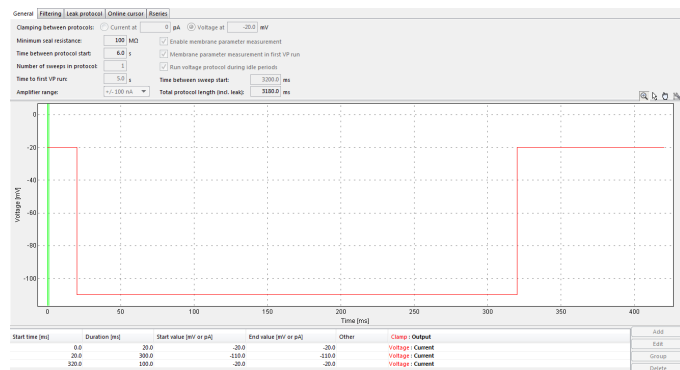
#1 Step IV protocol



#2 Ramp IV protocol



#3 Short pulse protocol



References:

1. Hsueh-Kai Chang, Jay-Ron Lee, Tai-An Liu, Ching-Shu Suen, Jorge Arreola, and Ru-Chi Shieh (2010). The Extracellular K⁺ Concentration Dependence of Outward Currents through Kir2.1 Channels Is Regulated by Extracellular Na⁺ and Ca²⁺. *J. Biol. Chem.* 285, 23115–23125.
2. Preisig-Muller R, Schlichthorl G, Goerge T, Heinen S, Bruggemann A, Rajan S, Derst C, Veh RW & Daut J (2002). Heteromerization of Kir2.x potassium channels contributes to the phenotype of Andersen's syndrome. *Proc Natl Acad Sci USA* 99, 7774–7779.

#4 Long ramp protocol

