

Application Report

Reliable $K_v2.1$ assays and pharmacology on the QPatch

$K_v2.1$ currents on the QPatch are stable and pharmacology values for the tested compounds are consistent with the literature.

Summary

- $K_v2.1$ is a major player in diabetes, neurodegeneration and other diseases
- Reliable currents on single-hole QPlates
- Consistent half-activation potentials for activation and inactivation
- Pharmacology values in agreement with literature values

Introduction

The voltage-gated potassium channel $K_v2.1$ is the main contributor to the delayed rectifier potassium current in neurons and in the cortex (Du et al. 1998, Murakoshi et al. 1999).

It has been suggested that $K_v2.1$ activation confers neuroprotection in a state of neuronal hyperexcitability via hyperpolarizing the membrane potential (Du et al. 2000, Misonou et al. 2005, Misonou et al. 2005a). Therefore, $K_v2.1$ may be an attractive target in epilepsy research. Also, the ion channel has previously been shown to play a central role in neurodegeneration, ageing and pancreatic β cell signaling.

Results and discussion

Activation of $K_v2.1$ results in stable currents

Cells were measured on single-hole QPlates in standard solutions (see Materials & Methods). The channel's current-voltage curve was determined by applying a voltage step protocol and measuring the resulting peak current (Fig.1 & 2).

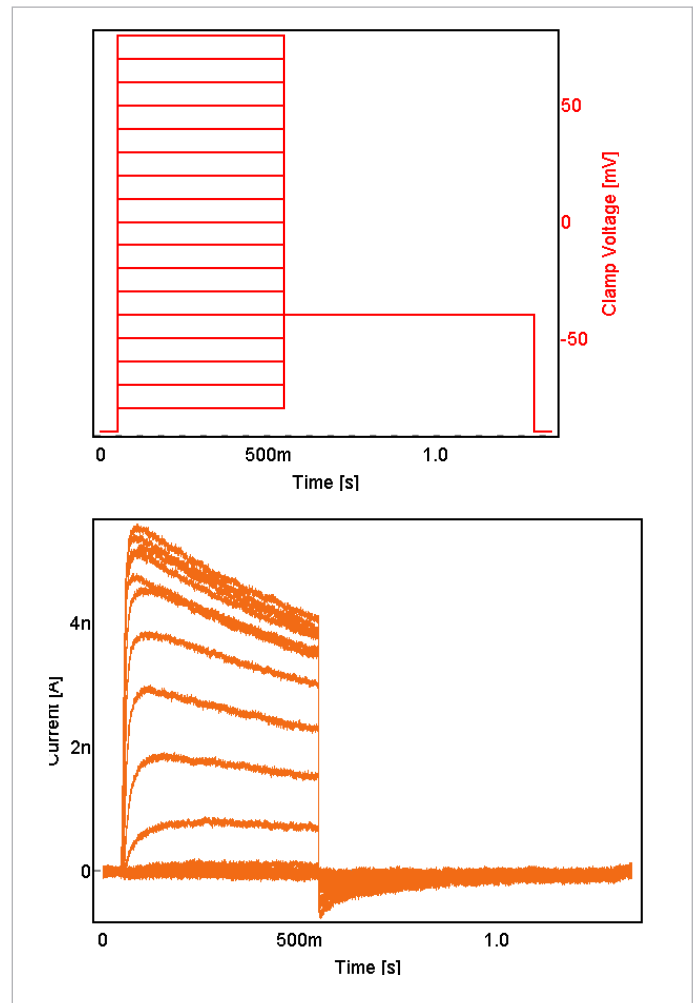


Fig. 1: CHO- $K_v2.1$ cells were depolarized from -90 mV to +80 mV in 10 mV steps (top). An example trace resulting from this protocol is shown in the lower figure.

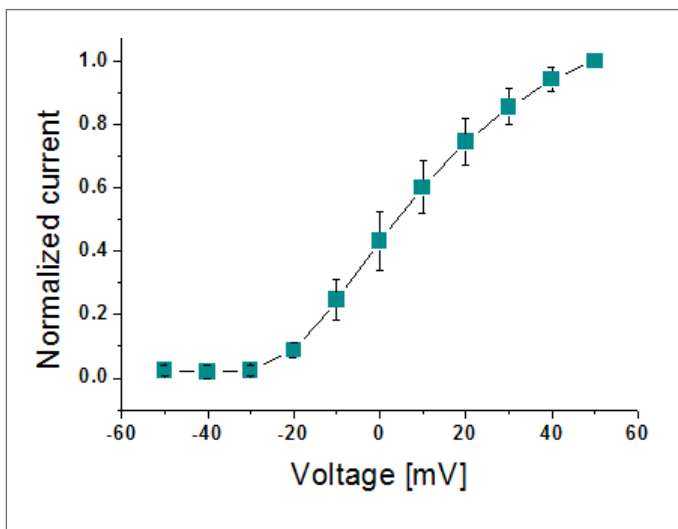


Fig. 2: Inactivation curve of $K_v2.1$. The current is normalized to the tail current amplitude at the -90 mV depolarization step. Data points represent normalized average tail current \pm SD, $n=18$.

Inactivation of $K_v2.1$

As before, CHO $K_v2.1$ cells were measured on single-hole QPlates in standard solutions (see Materials & Methods). A voltage step protocol was applied and a tail current was evoked by a subsequent depolarization step to +60 mV, where the resulting current was measured (Fig. 3 & 4). The average half-maximum inactivation potential ($V_{1/2inact}$) was -33 mV.

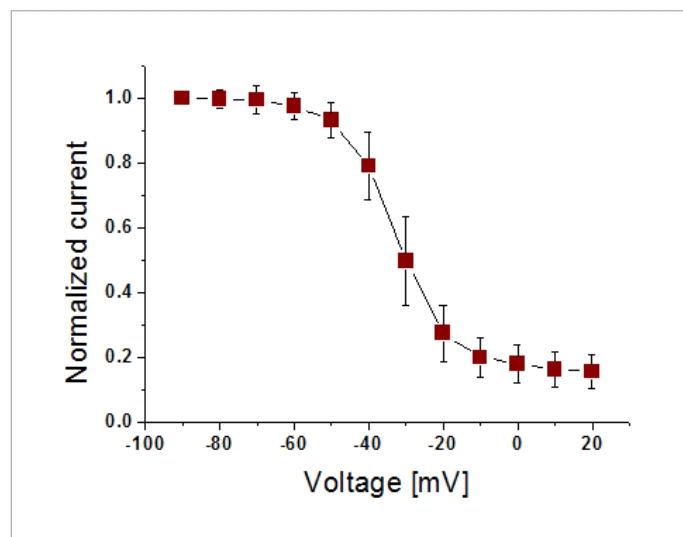


Fig. 4: Inactivation curve of $K_v2.1$. The current is normalized to the tail current amplitude at the -90 mV depolarization step. Data points represent normalized average tail current \pm SD, $n=18$.

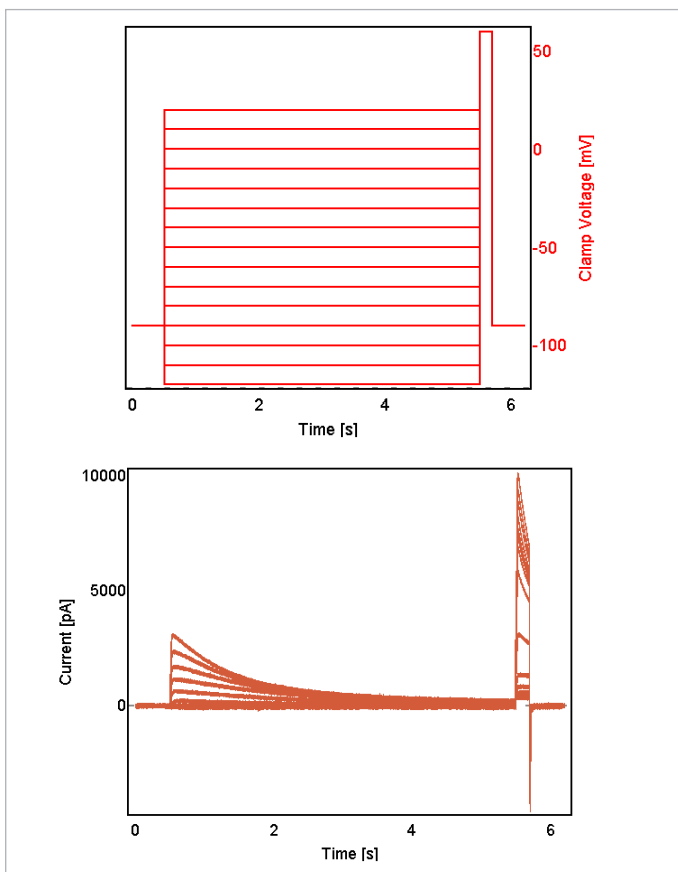


Fig. 3: CHO- $K_v2.1$ cells were clamped from -90 mV to +20 mV for 5 seconds in 10 mV increments, followed by a depolarization to +60 mV, at which the tail current was measured (top). An example trace resulting from this protocol is shown in the lower figure.

Pharmacology

CHO $K_v2.1$ cells were measured on single-hole QPlates in standard solutions (see Materials & Methods) with osmolarities adapted to the highest concentration of compound-containing solution (~350 mOsm). A depolarization from -80 mV to +20 mV was applied in saline solution, followed by cumulative addition of tetraethylammonium (TEA), nifedipine or saline. The lowest concentration of nifedipine was 1.2 μ M, followed by 3.7, 11.1, 33.3 and 100 μ M. The TEA solutions contained 0.37, 1.1, 3.3, 10 or 30 mM TEA. IC_{50} values of nifedipine (24 μ M) and TEA (17mM) (Fig. 5) are in agreement with literature values (Tagliatela et al. 1991, He et al 2006, Xian-Tao Li et al. 2015).

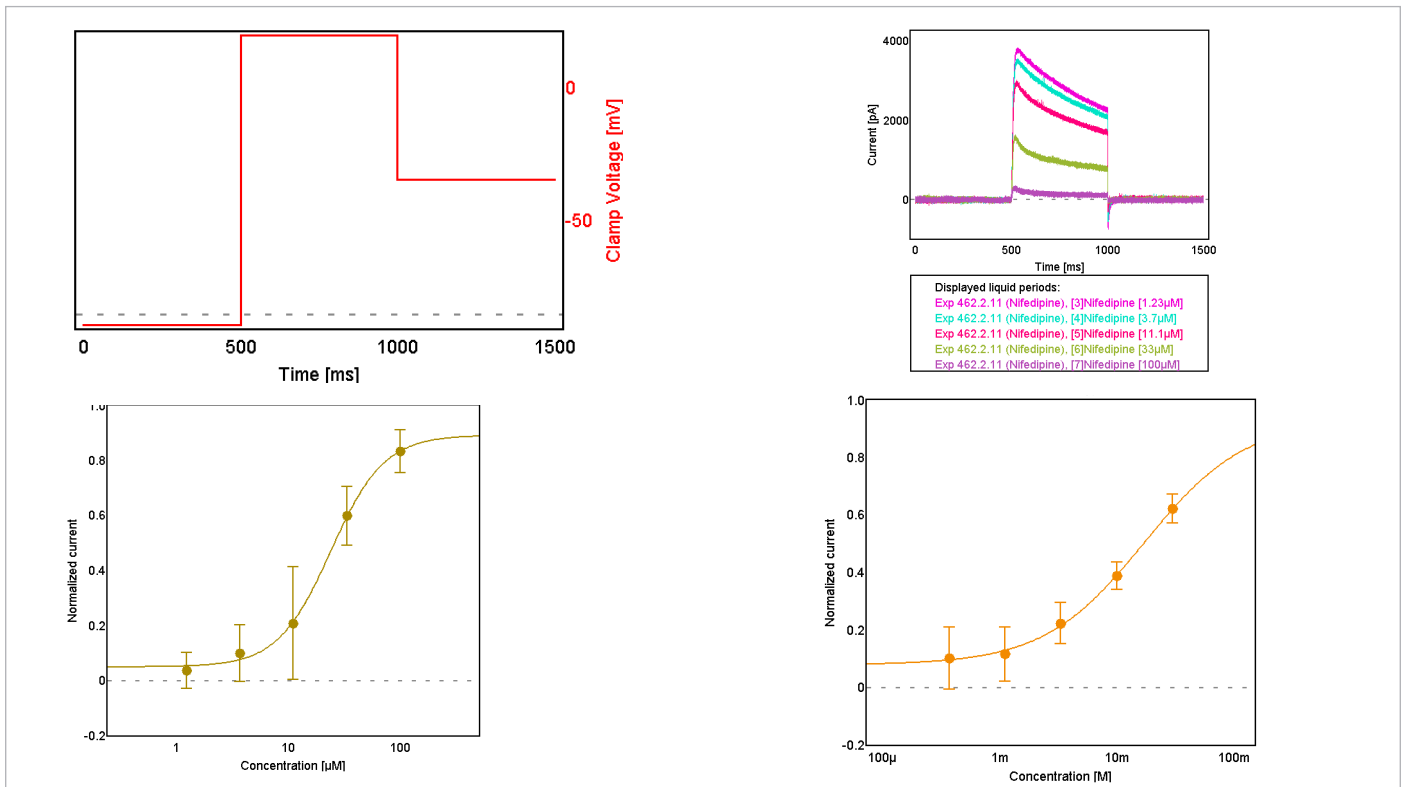


Fig. 5: Pharmacological inhibition of $K_v2.1$: Compound effects were assessed using a single depolarization step to +20 mV (left top). Both nifedipine (raw data example right top) and TEA inhibited $K_v2.1$ currents with IC_{50} values as displayed in the concentration-inhibition curve. A Hill equation was used to fit the data (solid lines). Data points represent normalized average current \pm SD, $n = 6$ (nifedipine) / 9 (TEA).

Methods

Cell culture for CHO cells, constitutively expressing human $K_v2.1$ (NP_004966.1), was performed according to B'SYS recommendations. Seeding and harvesting was done following Sophion's standard procedures.

For electrophysiological recordings, the extracellular solution contained (in mM): NaCl 145, HEPES 10, Glucose 10, KCl 4, $CaCl_2$ 2, $MgCl_2$ 2. pH=7.4 with NaOH, Osmolarity: 290 mOsm.

The intracellular solution was (in mM): Choline chloride 120, KOH/EGTA 31.25/10, HEPES 10, $CaCl_2$ 5.374, Na-ATP 4, $MgCl_2$ 1.75. pH=7.2 with KOH, Osmolarity: 280 mOsm.

All measurements were conducted with single-hole QPlates. The whole-cell was reached with a CHO standard protocol (1 second suction from -20 to -250 mbar with a subsequent pressure increment of -50 mbar until the whole-cell state is formed).

Conclusion

The depolarization of $K_v2.1$ cells, kindly provided by B'SYS, reliably resulted in potassium currents on the QPatch. The current-voltage relationship for activation, as well as inactivation

showed consistent values with minimal deviations. The QPatch is perfectly suited for the application of pharmacological compounds and the IC_{50} values for the drugs used in this assay align with literature values.

References:

- Du J, Tao-Cheng JH, Zerfas P et al., 1998. The K^+ channel, $K_v2.1$, is apposed to astrocytic processes and is associated with inhibitory postsynaptic membranes in hippocampal and cortical principal neurons and inhibitory interneurons. *Neuroscience*; 84(1):37-48.
- Du J, Haak LL, Phillips-Tansey E et al., 2000. Frequency-dependent regulation of rat hippocampal somato-dendritic excitability by the K^+ channel subunit $K_v2.1$. *J Physiol.*; 522 Pt 1:19-31.
- He Y, Kang Y, Leung YM et al., 2006. Modulation of $K_v2.1$ channel gating and TEA sensitivity by distinct domains of SNAP-25. *Biochem J.*; 396(2):363-9.
- Misonou H, Mohapatra DP, Menegola M et al., 2005. Calcium- and metabolic state-dependent modulation of the voltage-dependent $K_v2.1$ channel regulates neuronal excitability in response to ischemia. *J Neurosci.*; 25(48):11184-93.
- Misonou H, Mohapatra DP, Trimmer JS, 2005. $K_v2.1$: a voltage-gated K^+ channel critical to dynamic control of neuronal excitability. *Neurotoxicology*; 26(5):743-52.
- Murakoshi H, Trimmer JS, 1999. Identification of the $K_v2.1$ K^+ channel as a major component of the delayed rectifier K^+ current in rat hippocampal neurons. *J Neurosci.*; 19(5):1728-35.
- Tagliatalata M, Vandongen AM, Drewe JA et al., 1991. Patterns of internal and external tetraethylammonium block in four homologous K^+ channels. *Mol Pharmacol.*;40(2):299-307.
- Xian-Tao Li, Xiao-Qing Li, Xi-Mu Hu et al., 2015. The Inhibitory Effects of Ca^{2+} Channel Blocker Nifedipine on Rat $K_v2.1$ Potassium Channels. *PLoS One*; 10(4): e0124602.

Sophion Bioscience A/S, Baltorpevej 154, 2750 Ballerup, Denmark
Phone: +45 4460 8800 Fax: +45 4460 8899, E-mail: info@sophion.com

sophion.com