

# Cross-site study of neuronal K<sub>v</sub>7.2/7.3 heteromers on the Patchliner

Nanon Technologies GmbH, Munich and Apconix, Alderley Park, UK



## Summary

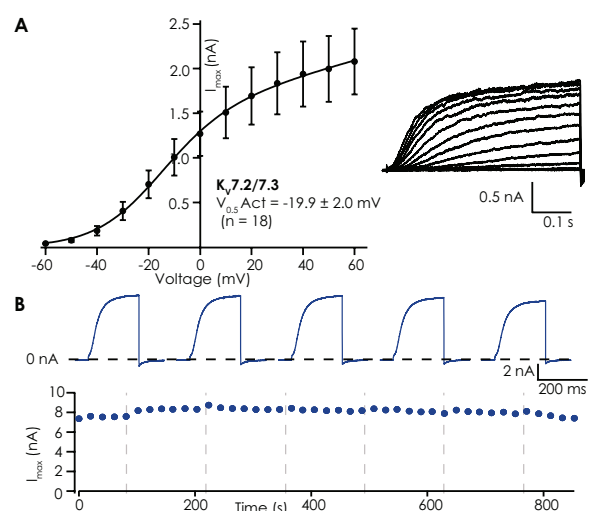
The K<sub>v</sub>7 channels are a family of voltage-gated potassium ion channels with five members (K<sub>v</sub>7.1 - 7.5) encoded by the KCNQ1-5 genes<sup>1</sup>. The channels exist as tetramers, with each subunit containing six transmembrane domains with cytoplasmic N- and C-termini. The long intracellular terminus is essential for tetramerization as well as interaction with critical regulators such as PIP<sub>2</sub>, calmodulin, protein kinase C and ankyrin G<sup>2</sup>. K<sub>v</sub>7-mediated currents are voltage activated, slowly activating and non-inactivating and are involved in repolarization of the cell membrane potential, thereby controlling cell excitability<sup>3</sup>. K<sub>v</sub>7.1 channels are primarily expressed in cardiac cells whereas K<sub>v</sub>7.2, K<sub>v</sub>7.3 and K<sub>v</sub>7.5 are widely distributed in neuronal and primary sensory cells<sup>2</sup>. K<sub>v</sub>7.2/7.3 heteromeric channels primarily underlie the neuronal M-current (I<sub>KM</sub>) which plays a crucial role in repolarizing neuronal membrane potential after a depolarizing input which limits repetitive firing and is, therefore, a key mechanism in spike frequency adaptation<sup>4</sup>.

Mutations in KCNQ2 and KCNQ3 have been shown to cause benign familial neonatal convulsions (BFNC), a rare inherited form of epilepsy, and neonatal epileptic encephalopathy (NEE)<sup>2,4</sup>. Mutations in the gene can affect voltage dependence of the resulting channel, alter the interaction with PIP<sub>2</sub>, affect the subcellular localization at key neuronal sites or cause loss of channel function<sup>2</sup>. Activation of K<sub>v</sub>7.2/7.3 is, therefore, a potential strategy to treat epilepsy with refigabine (ezogabine) approved as an adjunct therapy in partial-onset seizures in 2011, but was later withdrawn from

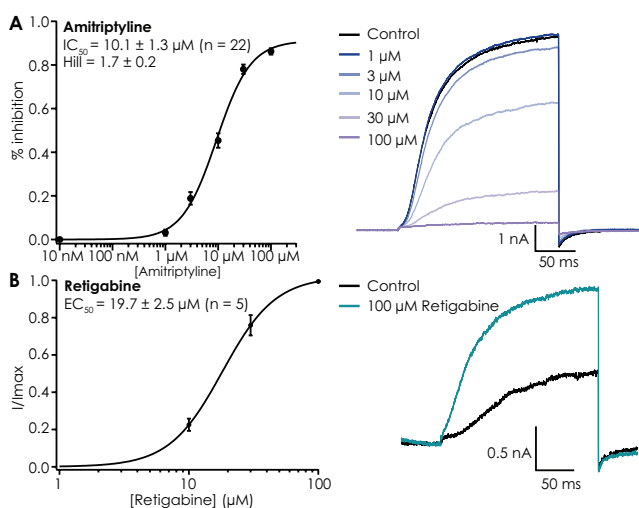
the market in 2017 due to adverse side effects. In addition to epilepsy, K<sub>v</sub>7.2/7.3 has also been proposed as a potential target to treat pain, neuropsychiatric disorders such as anxiety, attention deficit-hyperactivity (ADHD), mania, bipolar disorder and schizophrenia

## Results

K<sub>v</sub>7.2/7.3 stably expressed in HEK293 cells were used on the Patchliner at two different sites, Nanion HQ and Apconix. Current responses to increasing voltage steps were used to construct current-voltage plots. The curves were fit with a Boltzmann equation and the IV plot for an average of 18 cells recorded at Apconix is shown in Figure 1A with traces from an example cell shown in the inset. The average V<sub>half</sub> calculated from 18 individual plots was  $-19.9 \pm 2.0$  mV in excellent agreement with the literature<sup>6</sup>. At a second site (Nanon HQ), the average V<sub>half</sub> calculated from 16 individual plots was  $-22.2 \pm 1.8$  mV showing excellent consistency across the two sites. The recordings were also highly stable with little change in peak amplitude over a 15 minute measurement period. Figure 1B



**Figure 1:** **A** Current-voltage plot for K<sub>v</sub>7.2/7.3 with representative traces. **B** Stability of K<sub>v</sub>7.2/7.3 recordings. Currents could be recorded for over 15 minutes with little change in amplitude over time. Shown are raw traces from an exemplar cell and the corresponding time course in external solution.



**Figure 2: A** Amitriptyline inhibited  $K_{v7.2/7.3}$  in a concentration-dependent manner. The CRC for an average of 22 cells is shown and traces from an example cell. **B** Retigabine enhanced  $K_{v7.2/7.3}$ -mediated responses. The CRC for an average of 5 cells is shown and example traces from one cell +/- retigabine.

shows the timecourse and corresponding raw traces from an example cell.

Amitriptyline is a tricyclic antidepressant with analgesic and sedative properties, although it can be associated with severe side effects such as seizures<sup>7</sup>. It blocks both  $K_{v1.1}$  and  $K_{v7.2/7.3}$  channels which could underlie the neuroexcitatory side effects of the drug<sup>7</sup>. Amitriptyline blocked  $K_{v7.2/7.3}$ -mediated currents with an  $IC_{50}$  at one site (Nanion HQ; Figure 2A) of  $10.1 \pm 1.3 \mu\text{M}$  ( $n = 22$ ) and at the second site (Apconix) of  $6.9 \pm 1.2 \mu\text{M}$  ( $n = 8$ ) in excellent agreement with each other ( $P > 0.05$ , unpaired Student's  $t$  test) and with the literature<sup>7</sup>.

Retigabine is an anticonvulsant that enhances  $K_{v7.2/7.3}$ -mediated responses by shifting the voltage-dependence of activation to more hyperpolarized potentials<sup>9,10</sup>. In our experiments, retigabine enhanced  $K_{v7.2/7.3}$ -mediated responses in a concentration-dependent manner with an  $EC_{50} = 19.7 \pm 2.5 \mu\text{M}$  ( $n = 5$ ). Figure 2B shows the concentration response curve (CRC) for an average of 5 cells and the traces for an example cell in the presence and absence of  $100 \mu\text{M}$  retigabine.

In summary,  $K_{v7.2/7.3}$  expressed in HEK cells was recorded on the Patchliner at two different sites with very comparable results for  $V_{\text{half}}$  and pharmacology. The Patchliner is an excellent tool for studying  $K_{v7.2/7.3}$  and identifying novel anticonvulsants and analgesic compounds which enhance  $K_{v7.2/7.3}$  activity.

## References

1. Jepps, T.A., *et al.* 2021. *Front. Physiol.* 12:679317
2. Barrese, V., Scott, J.B., & Greenwood, I.A. 2018. *Annu. Rev. Pharmacol. Toxicol.* 58: 625 - 48

3. Tatulian, L., *et al.* 2001. 21 (15):5535 – 5545
4. Rogawski, M.A. *TINS.* 23(9): 393 - 398
5. Wulff, H., *et al.* 2009. *Nat. Rev. Drug Discov.* 8: 982 - 1001
6. Gamper, N., Stockand, J.D., & Shapiro, M.S. 2003. 23(1): 84 – 95
7. Punke, M.A., & Friederich, P. 2007. *Anesth. Analg.* 104(5): 1256 - 1264
8. Main, M.J. *et al.* 2000. *Mol. Pharmacol.* 58: 253 – 262
9. Wickenden, A.D., *et al.* 2000. *Mol. Pharmacol.* 58: 591 – 600

## Methods

### Cells

HEK293 cells stably expressing  $K_{v7.2/7.3}$  were used.

### Electrophysiology measurements

Cells were cultured and harvested according to Nanion's standard protocols. Cells were resuspended in external recording solution and stored in the CellHotel of the Patchliner before being dispensed into each well of the NPC-16 chip. Internal and external solution compositions are available upon request. For IV plots, voltage was increased in 10 mV steps from -60 mV to 60 mV from a holding potential of -80 mV. For amitriptyline, a single step to 60 mV from a holding potential of -80 mV was used and for retigabine, a single step to 0 mV was used. Data was analyzed in Igor or Prism. IV plots were fit with a Boltzmann equation and CRC plots with a Hill equation. Data is given as mean  $\pm$  S.E.M.

