

Structure-based identification of novel KNa1.1 inhibitors: a stratified target for *KCNT1*-related epilepsies



UNIVERSITY OF LEEDS

Bethan A. Cole¹, Rachel M. Johnson^{1,2}, Stephen P. Muench^{1,2}, & Jonathan D. Lippiat¹

¹School of Biomedical Sciences, ²Astbury Centre for Structural Molecular Biology, University of Leeds. UK.

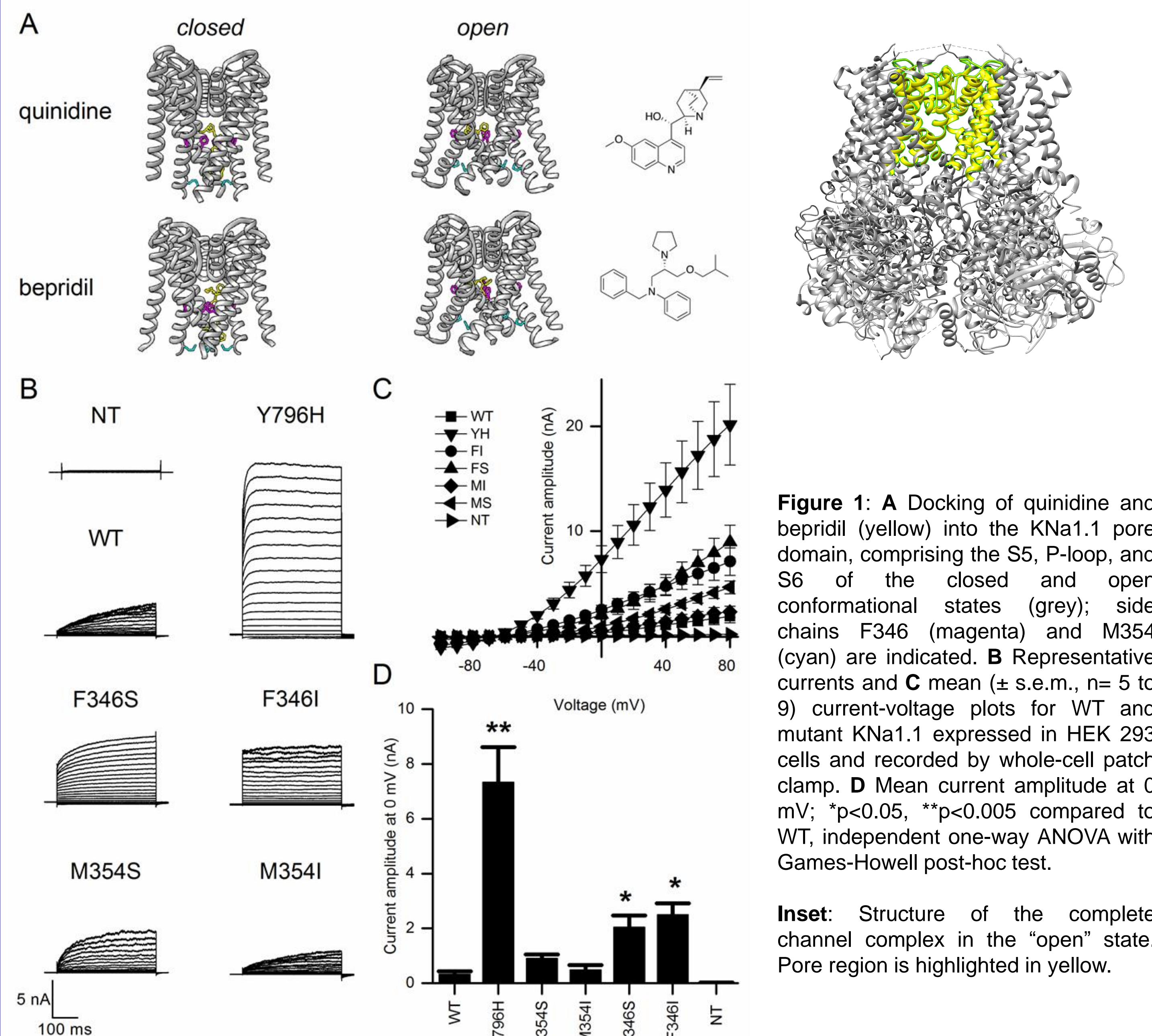
1. Background

- Missense, heterozygous, gain-of-function mutations of KNa1.1, a Na⁺-activated K⁺ channel encoded by *KCNT1* are associated with severe, pharmacoresistant epilepsies in children that are accompanied by psychomotor and intellectual disabilities. Malignant migrating partial seizures in infancy (MMPSI) and autosomal-dominant nocturnal frontal lobe epilepsy (ADNFLE) are two examples, but a number of other early-onset epileptic encephalopathies (EOEE) have also been identified [1,2,3]. There is a wide phenotypic spectrum associated with *KCNT1* mutations, and no specific inhibitors acting on the channel, making both prediction of disease outcome and treatment difficult [4]. The channel is widely distributed in the nervous system, and is thought to be involved in generation of the slow afterhyperpolarisation following a single action potential or trains of action potential firing [5].
- There are currently three known inhibitors of the channel: the antiarrhythmic drugs quinidine, bepridil, and clofilium [6,7]. All are non-selective and inhibit cardiac K⁺ and Na⁺ channels. Clinically, quinidine has had poor results, likely due to its non-selectivity and lack of potency. It is therefore important to develop a novel, selective inhibitor of KNa1.1.
- The structure of the chicken KNa1.1 channel was recently resolved using cryo-EM [8]. We hypothesised that this structure could be used to identify novel inhibitors *in silico* by docking a library of compounds into the pore-forming region of the channel. We have identified six novel compounds, all of which are more potent inhibitors of KNa1.1 than quinidine *in vitro*.

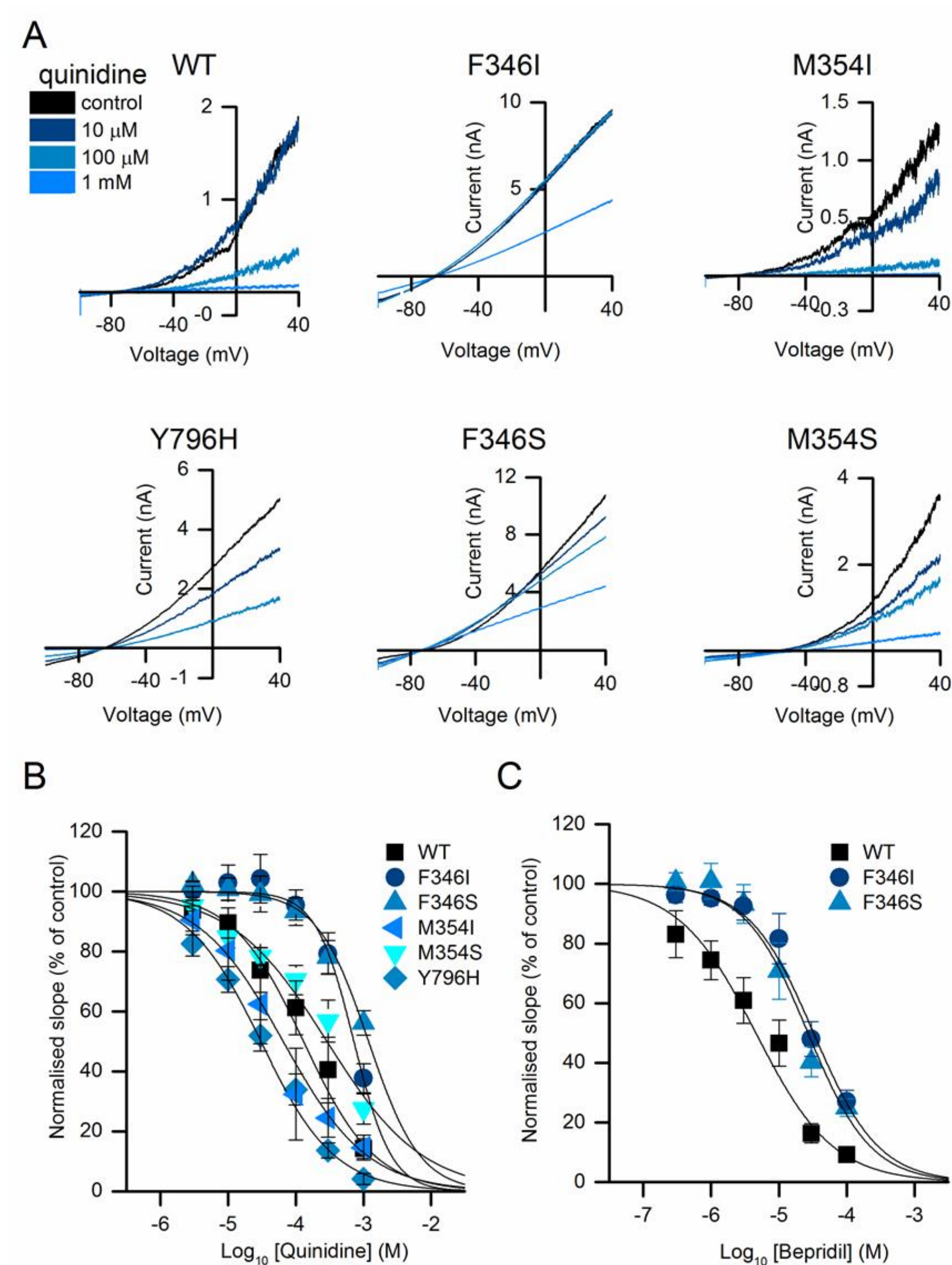
2. Methods

- Molecular docking was conducted using the cryo-EM structure of chicken KNa1.1, using SwissDock and GLIDE. Firstly, quinidine and bepridil were docked into the pore domain to identify potential binding sites, which were then used for virtual high-throughput screening using a Chembridge library of 100,000 commercially-available compounds. 17 high-scoring compounds were selected, ordered, and dissolved in DMSO (10 mM).
- HEK293-MSR cells were transiently transfected with wild-type (WT) or mutant human KNa1.1 constructs, or hERG.
- Currents were recorded by whole-cell patch clamp electrophysiology, using both voltage pulse and ramp protocols. Physiological solutions were used, with 10 mM Na⁺ included in the pipette (intracellular) solution for KNa1.1 recordings. Compound stock solutions were diluted with extracellular solution, and perfused onto cells serially or in increasing concentrations. Concentration-response curves were fitted using a Hill function, from which IC₅₀ values could be derived.
- Cell viability was assessed using a WST-1 assay of non-transfected (NT) HEK293 cells cultured in 96-well plates and measuring absorbance at 450 nm. Cell viability was determined using the equation:
% cell viability = ((A₄₅₀-A₆₅₀)_{experiment well} / ((A₄₅₀-A₆₅₀)_{control well}) x 100.

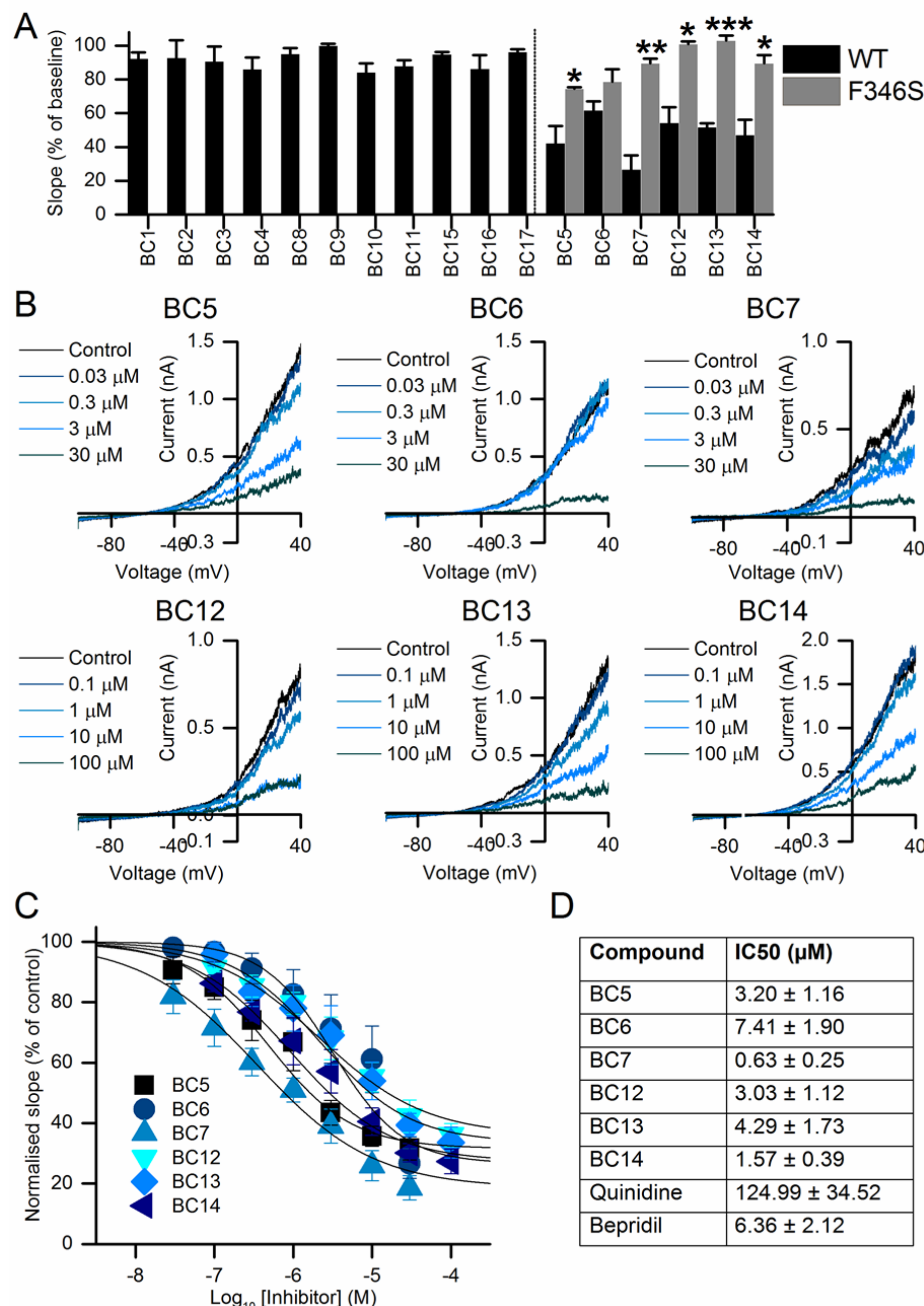
3. Results I



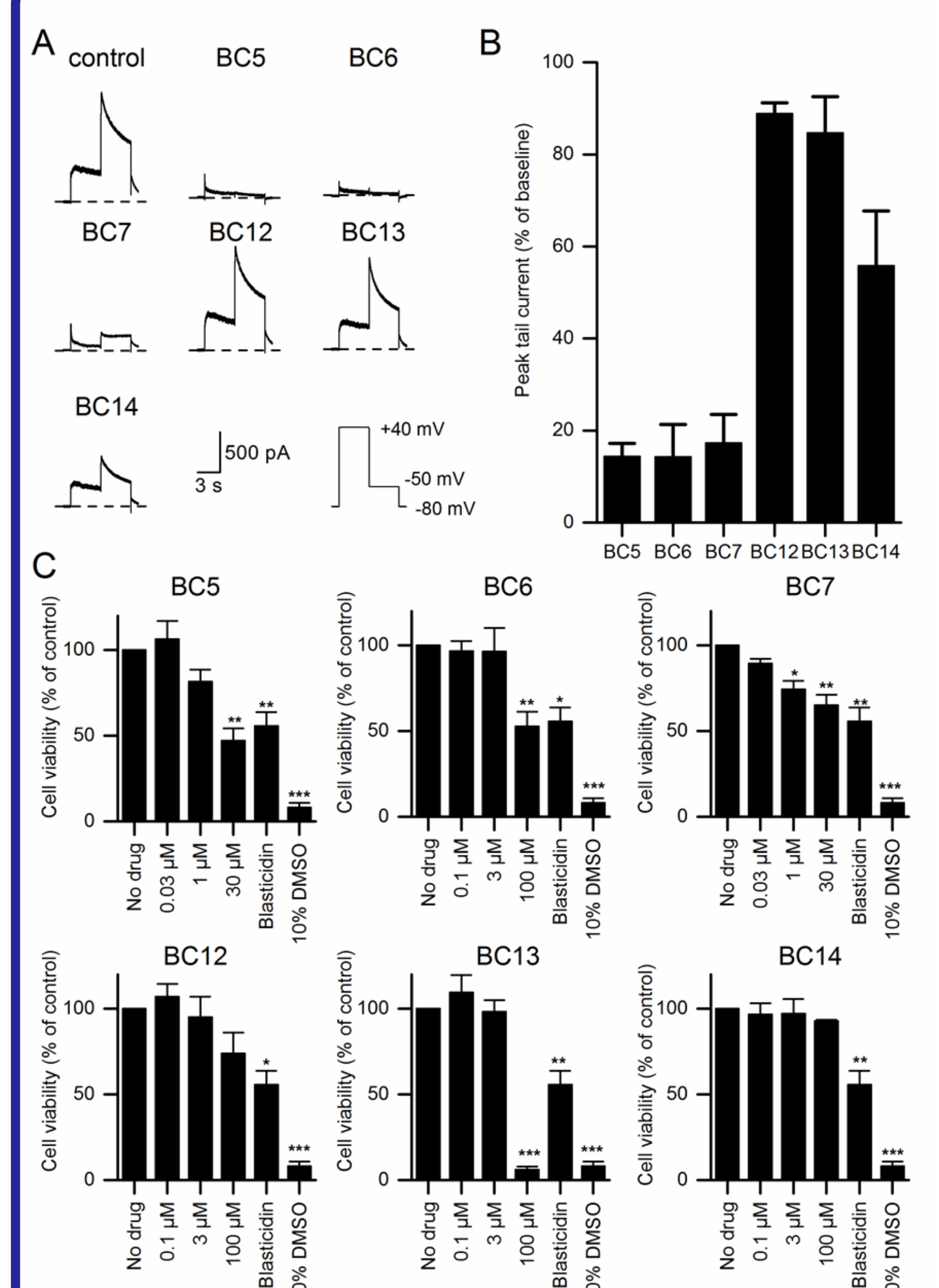
3. Results II



3. Results III



3. Results IV



4. Conclusions

- A phenylalanine residue, F346, in the pore-forming region of KNa1.1 is important for inhibition of the channel by quinidine and bepridil.
- Epilepsy-causing mutation Y796H increases quinidine sensitivity of the channel.
- The structure of the chicken KNa1.1 channel resolved by cryo-EM was successfully used to identify novel inhibitors of the channel using computer-aided methods.
- Reduced efficacy of the six compounds with F346S suggests they are specifically inhibiting the pore-forming region of KNa1.1.
- These are potential tool compounds or novel starting points for developing KNa1.1-specific inhibitors, though some may have toxic effects.

5. References/ acknowledgements

- Barcia et al (2012), Nat. Genet. 44 (11) 1255-9;
- Heron et al (2012), Nat. Genet. 44 (11) 1188-90;
- Lim et al (2016), J. Med. Genet. 53 (4), 217-25;
- Milligan et al (2014), Ann. Neurol. 75 (4) 581-90;
- Bhattacharjee et al (2003), J. Neurosci. 23 (37), 11681-91;
- Yang et al (2006) Neuropharmacology. 51 (4) 896-906;
- Rizzo et al (2016) Mol. Cell. Neurosci. 72, 54-63;
- Hite et al (2015) Nature. 527 (7577), 198-203.

We thank the BBSRC, Wellcome Trust, and Autifony Therapeutics for support.