

Co-expression of calcium and hERG potassium channels reduces the incidence of proarrhythmic events

Sara Ballouz ^{1,2,3,†}, Melissa M. Mangala ^{4,†}, Matthew D. Perry ^{2,4,†},
Stewart Heitmann ⁴, Jesse A. Gillis ³, Adam P. Hill ^{2,4,*}, and
Jamie I. Vandenberg ^{2,4,*}

¹Garvan-Weizmann Centre for Cellular Genomics, Garvan Institute of Medical Research, 384 Victoria Street, Darlinghurst NSW 2010, Australia; ²University of New South Wales, Sydney, Kensington, NSW 2052, Australia; ³Stanley Institute for Cognitive Genomics, Cold Spring Harbor Laboratory, Cold Spring Harbor, One Bungtown Road, NY 11724, USA; and ⁴Victor Chang Cardiac Research Institute, Lowy Packer Building, 405 Liverpool Street, Darlinghurst, New South Wales 2010, Australia

Received 16 June 2020; revised 25 August 2020; editorial decision 14 September 2020; accepted 17 September 2020

Aims

Cardiac electrical activity is extraordinarily robust. However, when it goes wrong it can have fatal consequences. Electrical activity in the heart is controlled by the carefully orchestrated activity of more than a dozen different ion conductances. While there is considerable variability in cardiac ion channel expression levels between individuals, studies in rodents have indicated that there are modules of ion channels whose expression co-vary. The aim of this study was to investigate whether meta-analytic co-expression analysis of large-scale gene expression datasets could identify modules of co-expressed cardiac ion channel genes in human hearts that are of functional importance.

Methods and results

Meta-analysis of 3653 public human RNA-seq datasets identified a strong correlation between expression of *CACNA1C* (L-type calcium current, I_{CaL}) and *KCNH2* (rapid delayed rectifier K^+ current, I_{Kr}), which was also observed in human adult heart tissue samples. *In silico* modelling suggested that co-expression of *CACNA1C* and *KCNH2* would limit the variability in action potential duration seen with variations in expression of ion channel genes and reduce susceptibility to early afterdepolarizations, a surrogate marker for proarrhythmia. We also found that levels of *KCNH2* and *CACNA1C* expression are correlated in human-induced pluripotent stem cell-derived cardiac myocytes and the levels of *CACNA1C* and *KCNH2* expression were inversely correlated with the magnitude of changes in repolarization duration following inhibition of I_{Kr} .

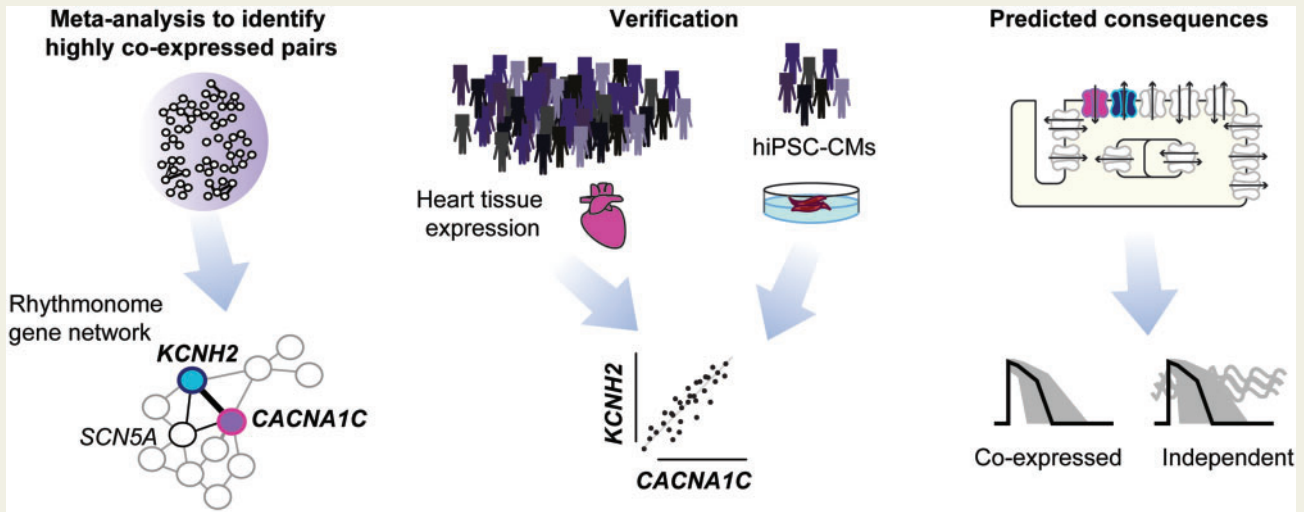
Conclusion

Meta-analytic approaches of multiple independent human gene expression datasets can be used to identify gene modules that are important for regulating heart function. Specifically, we have verified that there is co-expression of *CACNA1C* and *KCNH2* ion channel genes in human heart tissue, and *in silico* analyses suggest that *CACNA1C*–*KCNH2* co-expression increases the robustness of cardiac electrical activity.

*Corresponding authors. Tel: +61 2 9295 8771; fax: +61 2 9295 8601, E-mail: j.vandenberg@victorchang.edu.au (J.I.V.); E-mail: a.hill@victorchang.edu.au (A.P.H.)

[†]The first three authors contributed equally to the study.

Graphical Abstract



Keywords

Calcium channels • Potassium channels • Gene expression • Computational modelling • Cardiac arrhythmia

1. Introduction

Robust electrical signalling in the heart is critical for co-ordinating the efficient pumping of blood around the body. Failure of cardiac electrical signalling, even for just a few minutes, can have fatal consequences, with sudden cardiac death accounting for up to 10% of deaths in our community.¹ While we now have therapies for reducing the incidence of proarrhythmic triggers (e.g. beta-blockers²) and devices for terminating potentially fatal arrhythmias when they occur (e.g. implantable cardiac defibrillators³), predicting in advance who is more or less susceptible to sudden cardiac death and so who would most benefit from these therapies has proved to be extraordinarily difficult.⁴

The action potential (AP) of excitable cells, such as cardiac myocytes and neurons, reflects the orchestrated activity of at least a dozen distinct ion channels and electrogenic transporters. In such complex systems, with so many different molecular players, and hence degrees of freedom, it is not surprising that there are multiple combinations of different ion channel expression levels, that can repolarize the cell equally well.^{5,6} This has led to a paradigm shift in computational cardiology that relies not on generating idealized models based on mean data but rather development of populations of models that account for the observed variability in molecular inputs.^{7–9} A corollary of this concept is that while there are multiple possible ‘good-enough solutions’ that satisfy baseline requirements, not all combinations of ion channel expression will prove equally robust when faced with pathological insults.⁶

There is considerable heterogeneity of ion channel expression both within and between hearts.⁶ This heterogeneity, however, is not completely random. Rather there are reported examples of pairs or small modules of genes whose level of expression co-vary. For example, Bányász *et al.*¹⁰ and Rees *et al.*¹¹ demonstrated that the expression of I_{CaL} is correlated with the sum of the major repolarizing ion currents in guinea-pig and mouse, respectively. Furthermore, Rees and colleagues showed that in mice, this module of gene expression was important for

reducing the variability in the magnitude of the calcium transient. To date, there have been no similar studies in human heart tissue. Furthermore, the human heart expresses a different complement of repolarization channels compared to rodent hearts.¹² So even if this pattern was conserved, it is likely that different potassium channels would be involved in the human heart.

Outside the cardiovascular field, there is a considerable literature looking at the identification of co-expression networks among large-scale gene expression datasets.^{13,14} Such networks have been shown to encode functional information,¹⁵ with co-expression reflecting coregulation and co-functionality.^{11,16} Furthermore, meta-analytical methods have been developed, which aggregate large numbers of individual networks across multiple independent experiments to average away noise and remove bias that may be inherent in any one experimental approach thereby reinforcing those correlations that reflect real signals.^{17,18}

The aim of this study was to test the hypothesis that meta-analytical co-expression analysis of large-scale gene expression datasets would identify modules of co-expressed cardiac ion channel genes that are of functional importance for the human heart. This approach identified multiple pairs of co-expressed genes. To investigate whether pairs of co-expressed genes identified could be functionally significant, we focused on one co-expressed pair: *CACNA1C* (encodes the L-type calcium current, I_{CaL}) and *KCNH2* [encodes the rapid component of the delayed rectifier potassium current (I_{Kr}) which is often referred to as the human ether a-go-go related gene]. Using *in silico* approaches to generate populations of models of cardiac electrical activity with and without constrained patterns of co-expressed genes, we show that co-variation of I_{CaL} and I_{Kr} reduced the emergence of proarrhythmic early afterdepolarizations (EADs) and this protection persisted in the face of highly variable expression of other ion channels. Our modelling data also suggested that hearts with high expression of I_{CaL} and I_{Kr} are likely to be more susceptible to EADs when exposed to drugs that block I_{Kr} , the underlying basis

of drug-induced long QT syndrome,¹⁹ whereas low expression of I_{CaL} and I_{Kr} is associated with the greatest prolongation of repolarization duration without generation of EADs. This later prediction was verified in human cardiac myocytes derived from induced pluripotent stem cells. Together, our data, confirm that meta-analytic approaches of large-scale gene expression datasets are a valuable means for identification of gene modules that are important for regulating heart function.

2. Methods

A detailed version of the materials and methods section appears in the [Supplementary material online](#) that accompanies this article.

2.1 Analysis of public RNA-seq datasets

An aggregate co-expression gene network was built from public data, similar to that described previously.¹⁸ Briefly, 75 human RNA-seq expression experiments (listed in [Supplementary material online, Table S1](#)) that passed quality control and had a least 10 samples (3653 samples in total) were downloaded from the Gemma database.²⁰ (see [Supplementary material online](#) for more details). A co-expression network was generated for each experiment by calculating Spearman's correlation coefficients between every gene pair and then ranking these values.¹⁸ An aggregate gene co-expression network was then generated by averaging across all the individual networks, and re-ranking the final network. This final aggregate network was then used to determine the co-expression ranking between genes that encode for the set of ion channels and calcium handling proteins that determine the shape and duration of the human ventricular AP, the so-called rhythmome gene subset ([Supplementary material online, Table S2](#)).

2.2 GTEx dataset

Gene expression data from v8 of the GTEx atlas²¹ were downloaded from their webportal (<https://gtexportal.org/home/>). The dataset is human post-mortem RNA-seq across 31 tissues from 948 individuals, and included 429 atria samples and 432 samples from the left ventricle. Given that there may be significant mRNA degradation during the time between death and harvesting human heart samples, we repeated our analysis using only those samples with RNA integrity (RIN) scores >7.0; which corresponded to 472 heart samples (237 atria, 235 left ventricle).

2.3 Human-induced pluripotent stem cell-derived cardiac myocytes

Human iPSC lines, generated from healthy patients by Stanford Cardiovascular Institute Biobank, as previously described,²² were a generous donation from Joseph Wu (Stanford Cardiovascular Institute). HiPSC colonies were maintained on Matrigel[®] (Corning) coated plates in chemically defined medium (mTeSR1[™], StemCell technologies), and passaged using Dispase (StemCell technologies). HiPSC were differentiated using protocols as described previously (Perry 2020).²³ The CMs were seeded on a Matrigel-coated 96-well plate (Greiner Bio-One) and maintained in CM maintenance medium for 10–15 days before they were harvested for mRNA expression analysis. Total RNA was obtained from 40 000 hiPSC-CMs lysed using QIAzol Lysis reagent (Qiagen). RNA was purified using miRNeasy[®] Mini Kits (Qiagen) and analysed using Agilent Bioanalyzer pico-chip. RNA samples with RIN values >7.5 were hybridized with probes designed to detect 35 known rhythmome genes using nCounter (NanoString Technologies), which was performed at the Ramaciotti Centre for Genomics (UNSW). For

quantitative comparisons of different mRNA species across different runs, we normalized mRNA levels, relative to GAPDH and HPRT1, to account for any variation in mRNA loading levels.

Extracellular field potentials were recorded from hiPSC-CM monolayers using a Maestro-APEX multi-electrode array (MEA) system (Axion Biosystems, Atlanta, GA, USA) as previously described.²⁴ Field potential durations (FPD) were measured from the peak depolarization spike to the peak of the repolarization wave. After 30 min of pacing equilibration, electrograms were recorded for 5 min in control conditions, 5 min in the presence of cisapride (200 nM), and for 5 min in the presence of cisapride (200 nM)+nifedipine (50 nM). Measurements were averaged over the last 10 beats of each recording period. Twelve wells were plated for each line and all wells from which signals were recorded ($n = 7–12$) were included in the analysis. After the MEA recordings were completed, cells were harvested for mRNA analysis, as described above.

2.4 Statistical analysis

Summary data are shown as box and whiskers (median, box: 25–75%, extremities: 5–95%). Correlation coefficients for linear regressions were calculated in Prism (GraphPad Software, San Diego, CA) and statistical significance for whether the slope of the line was significantly different to zero was determined using the *F*-test.

2.5 Computer modelling

Human cardiac APs were simulated using a modified version of the endocardial configuration of the O'Hara-Rudy (ORD11) model,²⁵ with key conductances modified so that the model could reproduce the clinical phenotypes of long QT syndrome Types 1, 2, and 3 and maintain normal calcium transients as described by Mann *et al.*²⁶ and Krogh-Madsen *et al.*²⁷ The code was adapted to run in the Brain Dynamics Toolbox for Matlab.²⁸ To incorporate population variability in ion channel expression levels the maximum conductance for each current was multiplied by conductance scalar (G_x), that was drawn from a random log-normal distribution,²⁹ with unit mean and variance that was systematically manipulated from 0.05 to 0.5. All models were paced at 1 Hz with a stimulus of -40 mV and duration of 1 ms and allowed to equilibrate for 300 beats. We then analysed the next four beats (to allow for the possibility of development of alternans) after the equilibration stage. Individual beats were classified as ectopic if they had secondary peaks that were separated by >100 ms. The set of four beats were further classified as *alternans* if the profile of any of the APs deviated from each other by >1 mV at any time point. In a second set of simulations, we repeated the same method as above except that the random multipliers applied to both I_{CaL} and I_{Kr} were identical. This case, we denote co-expression, whereas the former case, we denote independent expression.

3. Results

3.1 Meta-analysis of cardiac ion channel gene expression

The shape and duration of APs in cardiac myocytes are determined by the orchestrated activity of voltage-gated sodium, calcium, and potassium channels, as well as a series of electrogenic transporters that regulate intracellular ion concentrations. These channels, transporters, and related intracellular calcium handling proteins are encoded by a few dozen genes, sometimes referred to as the rhythmome³⁰ (a full list is provided in [Supplementary material online, Table S2](#)).

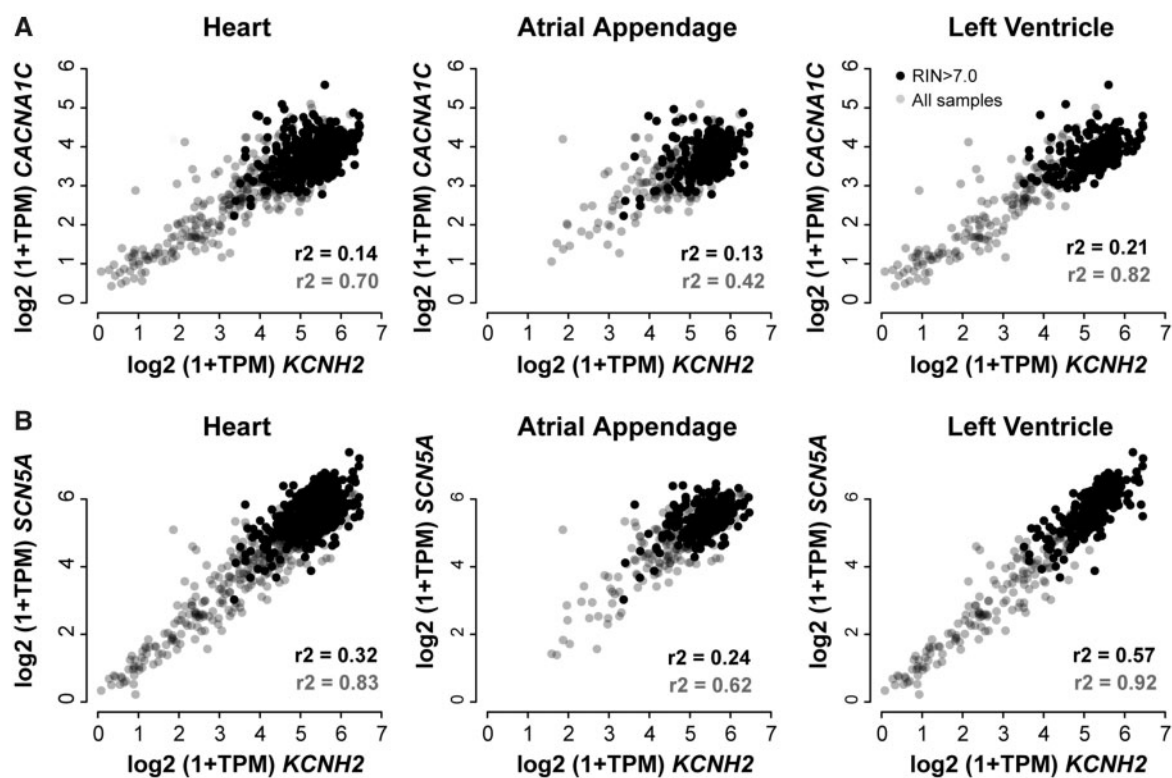


Figure 2 Expression levels of (A) *CACNA1C* and (B) *SCN5A* vs. expression levels of *KCNH2* in the GTEx dataset. Expression levels are expressed as $\log_2(\text{transcripts per million} + 1)$. Data points for samples with a RNA integrity number (RIN) >7.0 are shown in black (237 atria, 235 left ventricle samples) and all other sample data points are shown in grey (192 atria, 197 left ventricle samples). The left panel shows all heart samples, the middle panel shows atrial appendage samples, and the right panel shows left ventricular samples. Correlation coefficients shown in black and grey at the bottom right of each panel correspond to the full dataset (grey) and the samples with a high RIN (black). Correlation coefficients for linear fits to the data are shown in each panel. It is notable that correlations are stronger for the ventricular samples than for atrial samples.

to be important contributors to repolarization in the human heart.³² Third, previous studies in rodents have shown that there is co-expression of I_{CaL} and repolarization currents,^{10,11} suggesting that this may be a relationship conserved across species. And finally, there are highly specific pharmacological tools to enable specific dissection of the functional role of these channels.³³

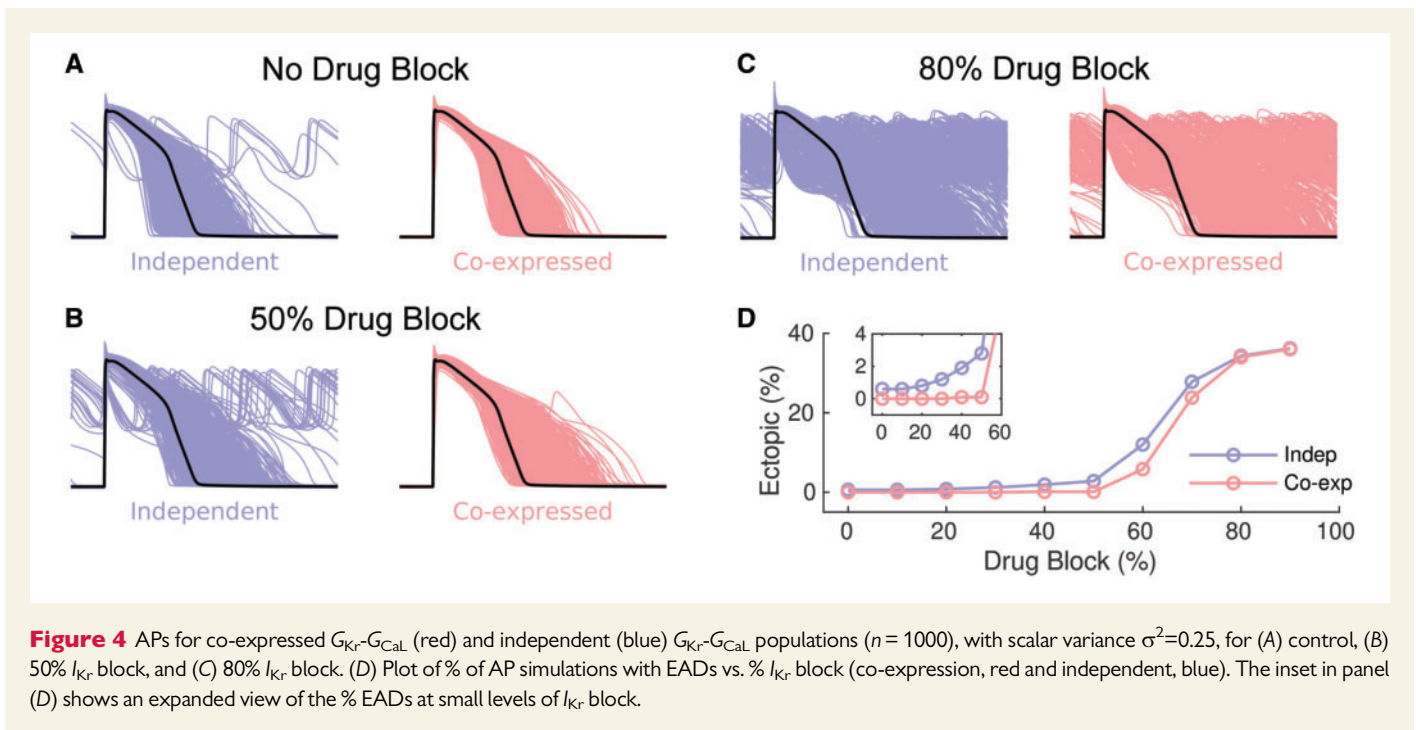
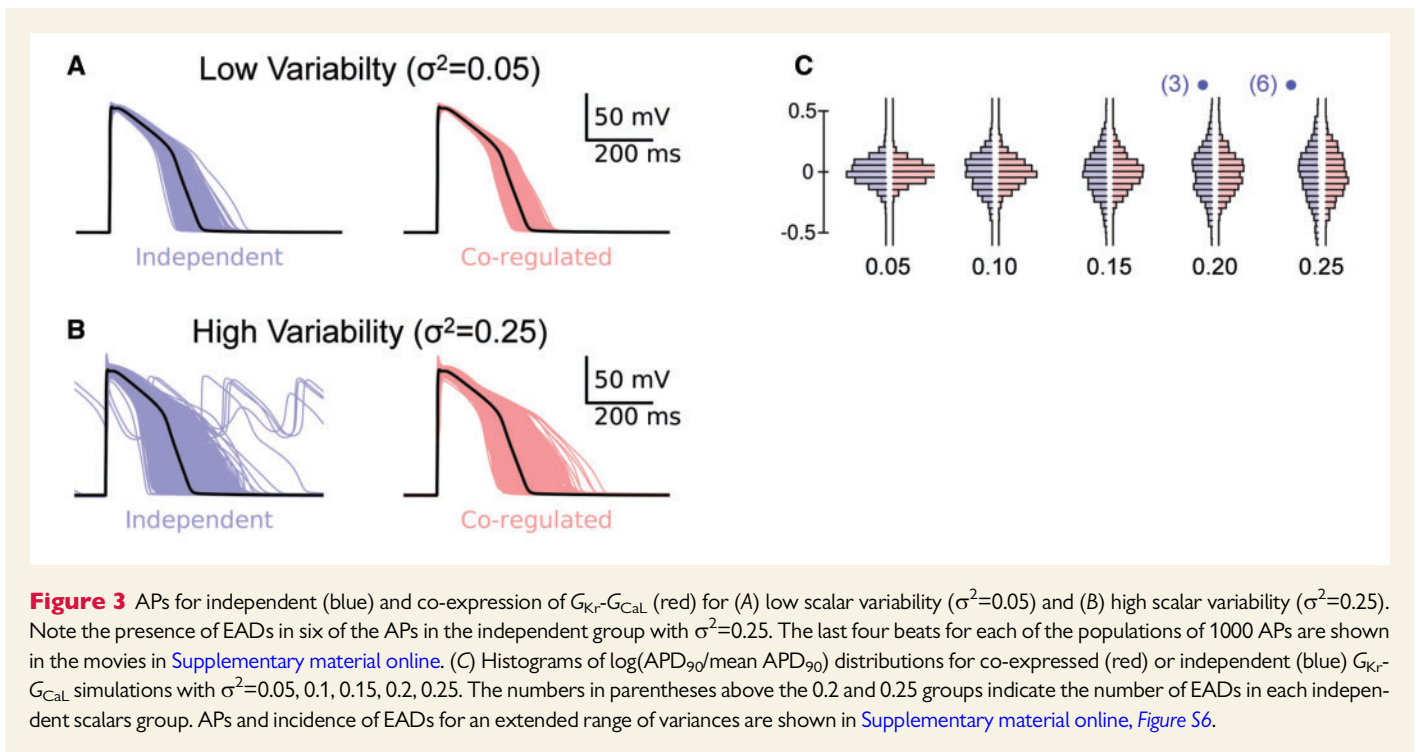
3.2 Co-expression of *CACNA1C* and *KCNH2* in human heart tissue

The public RNA-seq datasets included in our analyses spanned a wide range of tissue types and different experiments. This was a deliberate strategy to minimize the confounder effects that may be seen in any one type of experimental approach. Unfortunately, there are too few human heart samples to undertake specific co-expression meta-analysis just on heart samples. So, to investigate whether these co-expression patterns were observed in human heart, we looked at the GTEx dataset of post-mortem tissue samples.²¹ While we acknowledge there are limitations associated with the GTEx dataset, for example, the majority of samples is from diseased hearts and there is a significant variation in the time between death and tissue collection, it is clear that there is a correlation between the expression of *CACNA1C* and *KCNH2* (as well as between

SCN5A and *KCNH2*) (Figure 2). Furthermore, this correlation was much stronger in the ventricular samples than in the atrial samples. It is also worth noting that in other tissues, collected as part of the GTEx dataset, that show robust expression of both *KCNH2* and *CACNA1C*, the correlation between expression of *KCNH2* and *CACNA1C* was also observed (e.g. oesophagus and colon; see Supplementary material online, Figure S3).

3.3 *In silico* analysis of impact of co-expression of *CACNA1C* and *KCNH2* on cardiac electrical activity

Given the correlation between *CACNA1C* and *KCNH2* expression was most apparent in ventricular tissue, and these two genes encode for the most important currents (I_{CaL} , inward current and I_{Kr} , outward current) controlling repolarization in the human ventricular AP under resting conditions,²⁹ we next investigated how the co-expression of these two genes would influence ventricular repolarization. Initially, we used an *in silico* approach based on the modified O-Hara-Rudy model of the human ventricular myocyte with the baseline conductance scalars modified so that the model could reproduce the clinical phenotypes of long QT



syndrome Types 1, 2, and 3 and maintain normal calcium transients as previously described^{26,27} ([Supplementary material online, Figure S4](#)).

First, we simulated a population of 1000 human cardiac APs where the ion conductances for each cell were scaled randomly using a log-normal distribution, that is, meaning the probability of the conductance being doubled was the same as that of it being halved (see [Supplementary material online, Figure S5](#) for details of generation of

scalars). The population of models generated when all conductance scalars were varied independently ([Supplementary material online, Movie S1](#)) are denoted independent in [Figure 3](#). We then repeated the simulation but forced the conductance scaling factors for I_{K_r} and I_{CaL} to be identical to each other, while the scaling factors for all other conductances were allowed to vary independently, denoted co-expression in [Figure 3](#) (and see [Supplementary material online, Movie S2](#)).

which is the predominant cause of drug-induced long QT syndrome.¹⁹ Populations of APs obtained for independent and co-expression populations with I_{K_r} block of 0%, 50%, and 80% are illustrated in Figure 4. As the extent of I_{K_r} block is increased (from A to C), the proportion of APs producing EADs increases. It is also clear that at lower levels of I_{K_r} block, EADs were more frequent when G_{CaL} and G_{K_r} scalars were modulated independently (see Figure 4B and inset to Figure 4D). However, the proportion of simulations developing EADs in both the independent and co-expression populations becomes similar when the extent of I_{K_r} block exceeds 80% (Figure 4D).

It is well established that the extent of repolarization prolongation seen in patients prescribed drugs that block I_{K_r} is highly variable and only a subset of patients exposed to drugs that block I_{K_r} ³⁴ or with a mutation causing 50% loss of I_{K_r} function,³² will develop life-threatening arrhythmias. This is consistent with the prediction made by our simulated drug block experiments shown in Figure 4. We therefore asked whether the data from the co-expression datasets could tell us anything about what factors might predispose to lesser or greater extent of AP prolongation and/or the development of EADs in the presence of a drug that blocks I_{K_r} . Analysis of the subset of scalars within the co-expression dataset that produced the 50 longest APDs without EADs, compared to the subset of scalars that resulted in APs with EADs, is illustrated in Figure 5A and B, respectively. Notably, the longest APs without EADs had low G_{CaL} scalars (and hence low G_{K_r} scalars before addition of drug block). Conversely, the APs that developed EADs had higher G_{CaL} and G_{K_r} scalars. In Figure 5C, we have plotted the APD₉₀ values for cells in the highest (red) and lowest (blue) quartiles of G_{CaL} – G_{K_r} scalars. As expected, the low G_{CaL} group showed longer APD₉₀ values on an average compared to the high G_{CaL} group (see the continuous lines in Figure 5C). Furthermore, for the 70% I_{K_r} block scenario, 44% of the high G_{CaL} group have developed EADs whereas only 7% of the low G_{CaL} group have developed EADs (compare red and blue bars at 70% drug block in Figure 5C). Thus, higher G_{CaL} is associated with a greater risk of developing EADs in response to moderate levels of I_{K_r} block. A similar pattern of results was observed when G_{CaL} and G_{K_r} were allowed to vary independently, except that in this scenario the difference between the high G_{CaL} and low G_{CaL} groups was even more dramatic at lower levels of I_{K_r} block (Supplementary material online, Figure S7). A corollary of our prediction that patients with high G_{CaL} are more susceptible to EADs when exposed to a drug that blocks I_{K_r} , is that co-administration of a drug that blocks I_{CaL} would reduce the incidence of EADs. A drug that inhibited I_{CaL} by 20% caused a modest decrease in the percentage of cells with EADs and reduction of I_{CaL} by 50% had a correspondingly larger effect (Supplementary material online, Figure S8).

3.5 Validation of *in silico* predictions in human cardiac myocytes derived from induced pluripotent stem cells

So far, we have identified modules of co-expressed genes in RNA-seq datasets, demonstrated using *in silico* modelling of the human ventricular AP that co-expression of *CACNA1C* and *KCNH2* would produce more robust electrical signalling, and shown that the extent of prolongation of repolarization seen following I_{K_r} drug block should be inversely proportional to the level of *CACNA1C*/*KCNH2* expression. We next wished to investigate whether this could be verified in heart tissue, given that *KCNH2*-encoded I_{K_r} is the major repolarization current in human heart but not in rodents.¹² Due to the difficulties in obtaining adult human heart cells, we chose to use human-induced pluripotent stem cell-derived cardiac myocyte cell lines to determine whether the predictions

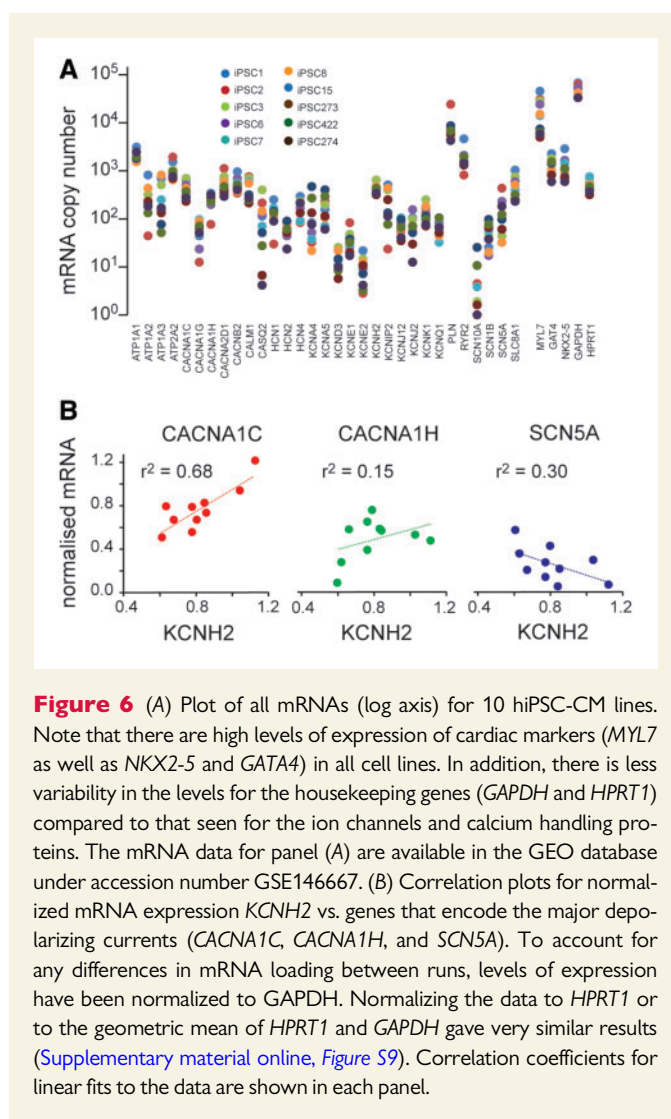


Figure 6 (A) Plot of all mRNAs (log axis) for 10 hiPSC-CM lines. Note that there are high levels of expression of cardiac markers (*MYL7* as well as *NKX2-5* and *GATA4*) in all cell lines. In addition, there is less variability in the levels for the housekeeping genes (*GAPDH* and *HPRT1*) compared to that seen for the ion channels and calcium handling proteins. The mRNA data for panel (A) are available in the GEO database under accession number GSE146667. (B) Correlation plots for normalized mRNA expression *KCNH2* vs. genes that encode the major depolarizing currents (*CACNA1C*, *CACNA1H*, and *SCN5A*). To account for any differences in mRNA loading between runs, levels of expression have been normalized to *GAPDH*. Normalizing the data to *HPRT1* or to the geometric mean of *HPRT1* and *GAPDH* gave very similar results (Supplementary material online, Figure S9). Correlation coefficients for linear fits to the data are shown in each panel.

from our *in silico* prediction holds up in human heart cells. First, we extracted mRNA from human-induced pluripotent stem cell-derived cardiomyocytes (hiPSC-CMs)³⁵ obtained from patients with no known heart disease. All samples contained high levels of cardiac marker genes (*MYL7*, *GATA4*, and *NKX2.5*) (Figure 6A). The levels of expression of most rhythmome genes showed variations between the samples that spanned approximately an order of magnitude. As these cell lines represent an embryonic phenotype, they have low levels of expression of a number of ion channel genes, including, for example, *SCN5A*, *KCND3*, and *KCNJ2*. However, they do express robust levels of *KCNH2* and *CACNA1C*. Most importantly, there was also a strong correlation between the level of expression of *KCNH2* and *CACNA1C* (Figure 6B, $r^2=0.68$).

We next measured the duration of repolarization from extracellular field potentials (FPD) recorded from monolayers of hiPSC-derived cardiac myocytes paced at 1 Hz and exposed to drugs that specifically block I_{K_r} , cisapride,³³ and I_{CaL} , nifedipine.³³ Cisapride (200 nM) caused a prolongation of repolarization that varied between 17.9 ± 7.4 ms (mean \pm SD, $n=12$) for the least sensitive cell line to 98 ± 41 ms (mean \pm SD, $n=11$) for the most sensitive cell line (Figure 7A). There was also a significant inverse correlation between the extent of FPD prolongation and the [*KCNH2*–mRNA] measured for each cell line (Figure 7B).

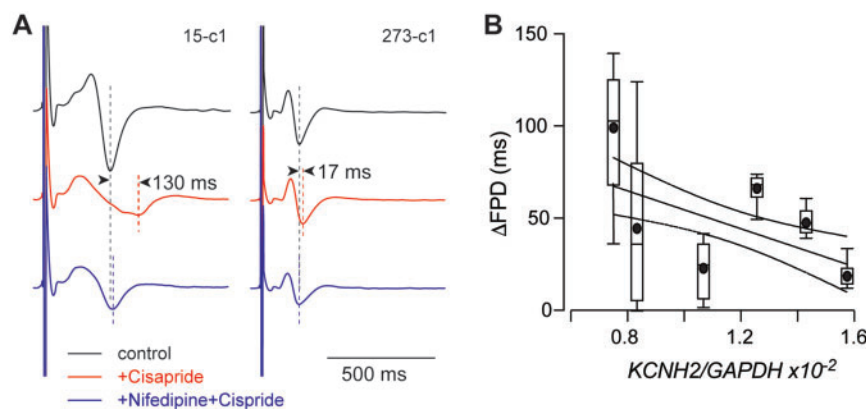


Figure 7 (A) Example extracellular electrograms recorded from two separate hiPSC-CM lines under control (black) conditions, after 5 min perfusion with cisapride (200 nM, red) followed by 5 min perfusion with cisapride (200 nM)+nifedipine (50 nM, blue). In cell line 15-c1, there was a 130 ms prolongation in field potential duration (FPD) which almost completely reversed on addition of nifedipine. In cell line 273-c1, there was a much smaller prolongation of 17 ms with addition of cisapride that again reversed on the addition of Nifedipine. (B) Summary of Δ FPD for six cell lines plotted against the level of *KCNH2* mRNA expression (normalized to *GAPDH* expression). Data shown as box and whiskers (median, box: 25–75%, extremities: 5–95%, mean: filled circle, $n = 7$ –12 replicates for each cell line). The line of best fit is shown as a solid line with 95% confidence intervals as dashed lines; the slope of the line is significantly different to zero (F -test, $P = 0.0010$). The micro-electrode array data for panels (A) and (B) are available in the DRYAD repository at doi: <https://doi.org/10.5061/dryad.612jm640k>. The mRNA data for panels (B) and (C) are available in the GEO database under accession number GSE146667.

Furthermore, the extent of FPD prolongation caused by cisapride in each cell line was almost exactly reversed by Nifedipine (50 nM) (Figure 7A).

4. Discussion

A better understanding of the molecular and cellular level control of cardiac electrical activity could pave the way for improved detection of patients at risk of cardiac arrhythmias. Here, we report that the levels of *KCNH2* and *CACNA1C* mRNA expression, in generic tissue RNA-seq datasets (Figure 1) and the GTEx Atlas of heart tissue samples (Figure 2), vary considerably from person to person but that the levels of *KCNH2* and *CACNA1C* expression are tightly correlated. Our *in silico* population modelling of the human ventricular AP shows that co-expression of I_{Kr} (encoded by *KCNH2*) and I_{CaL} (encoded by *CACNA1C*), in the context of random variations in the densities of all cardiac ion channels, results in a narrower range of APD₉₀ values at baseline (Figure 3) and helps to protect against the development of EADs (Figures 3 and 4). Our modelling approach also suggests that individuals with higher I_{CaL} levels will be more susceptible to EADs when exposed to drugs that block I_{Kr} (Figure 5), whereas the subset of individuals with lower levels of I_{CaL} will develop more prolongation of repolarization without developing EADs. Importantly, independent data obtained from human iPSC-derived cardiac myocytes have validated key predictions from our *in silico* analyses (Figure 7).

Over the last few years, numerous groups have shown that there is considerable heterogeneity of ion channel expression among excitable cells, including neurons⁵ and cardiac myocytes.^{6,10,36} Furthermore, since the pioneering work of Marder and colleagues, the presence of modules of co-expressed ion channel genes has also been well appreciated.¹⁴ These previous studies, however, relied on patch clamp analysis of isolated cells^{10,14} and qPCR analysis¹⁴ or single-molecule fluorescence *in*

situ hybridization³⁶ of individual ion channel mRNAs, which are highly laborious techniques and so have been restricted to only a few important ion channel genes. The advent of high-throughput transcriptomic analyses has greatly facilitated the identification of conserved networks of co-expressed gene modules among the entire set of expressed genes.¹⁸ By applying meta-analytic approaches to a large number of independent datasets one can more readily discern genuine co-expression signals from noise, as well as explore smaller gene modules.¹⁸ In our analysis of a subset of rhythmome genes within a large library of public RNA-seq datasets, we identified multiple correlations within the subset of genes that predominantly affects calcium handling, the subset that predominantly affects membrane potential as well as between these two subsets of genes (see Figure 1). The nodes within each of these clusters that are most closely connected within the rhythmome relative to all other genes are *KCNH2* and *CACNA1C* (Supplementary material online, Figure S1). This is analogous to a network of network,³⁷ where the *KCNH2*–*CACNA1C* link provides an interconnection of the two sub-networks. Independent evidence to corroborate an important link between calcium handling and regulation of cardiac AP duration comes from the large genome-wide association studies of QT interval duration which identified SNPs in a number of calcium handling genes as well as *KCNH2* as being important determinants of QT interval in the population.³⁸

Due to the large number of associations, we were testing for in our network analyses, it is possible that some would occur by chance. It was therefore critical that we investigated whether any of the identified correlations could be verified in human heart tissue and whether they were of functional importance. That we were able to confirm the presence of important co-expression modules in an independent dataset, that is, GTEx²¹ heart tissue samples (see Figure 2 and Supplementary material online, Figure S2) as well as in hiPSC-CM (see Figure 6) provides important corroborative evidence to support the conclusion that co-expression patterns are likely to be real and have physiological relevance. The patterns of co-expressed genes we observed also show some

important similarities to previous studies in rodents. Bányász et al.¹⁰ and Rees et al.¹¹ have both demonstrated that the expression of I_{CaL} is correlated with the sum of the major repolarizing ion currents in guinea-pig and mouse, respectively. However, the molecular players involved in cardiac electrical activity in rodents are quite distinct to humans.¹² In humans, the most important determinant of repolarization duration at resting heart rates in the absence of adrenergic stimulation is I_{Kr} ,²⁹ whereas in guinea-pig I_{Ks} and I_{Kr} play equally important roles¹⁰ and in mice the fast component of the transient outward current ($I_{to,f}$) and the ultra-rapid delayed rectifier (I_{Kur}) are the major repolarization currents.¹¹ Our study, however, is the first to demonstrate the importance of co-expressed genes important for repolarization in human heart tissue.

Identifying modules of co-expressed genes are the first step in seeking to understand the logic of complex systems.⁶ The next challenge is understanding how such modules impact function in health and disease. Recently, Rees et al.¹¹ have demonstrated that modules of co-expressed depolarization and repolarization currents can help to ensure normal amplitude calcium transients in murine heart cells, a critical determinant of overall heart function. We have extended these studies to show that, co-expression of *KCNH2* (repolarization) and *CACNA1C* (depolarization) channels not only limits the extent of variability in duration of the ventricular AP (Figure 3) it also helps to protect the heart from early after depolarizations when they are exposed to drugs that block I_{Kr} (Figure 5).

It is worth noting that I_{Kr} is close to fully activated during the cardiac AP at resting heart rates.³⁹ Conversely, I_{CaL} is only partially activated under resting conditions and Ca^{2+} fluxes through *CACNA1C* channels can increase markedly under conditions of adrenergic stress or exercise.³⁹ An unopposed increase in I_{CaL} would likely be proarrhythmic. However, adrenergic stress also results in a marked increase in K^+ flux through the slow delayed rectifier K^+ channels encoded for by *KCNQ1* and *KCNE1*.³⁹ Our meta-analysis data (Figure 1) do not indicate a close correlation between *KCNQ1/KCNE1* and *CACNA1C*, suggesting that these channels are not co-regulated at a transcriptional level. However, given that acute responses to adrenergic stress are mediated by post-translational modifications, it is likely co-regulation of I_{CaL} and I_{Ks} occurs at a post-transcriptional level. Indeed, there is considerable evidence for co-regulation of cardiac ion channels occurring at post-transcriptional levels.^{36,40} For example, recent work from Robertson and colleagues demonstrated co-translation of *SCN5A* and *KCNH2* channels in human iPSC-derived cardiac myocytes.³⁶ There is also evidence to suggest that different ion channels traffic together to reach similar destinations at the cell surface, for example, the co-assembly of I_{Na} , I_{to} , and I_{K1} at intercalated discs^{40–42} has an important influence on intercellular conduction. Thus, any complete cardiac electrophysiology modelling approach will have to include not just co-regulation at the level of expression but also co-regulation occurring at the post-transcriptional level.

Our *in silico* studies, incorporating co-expression of *KCNH2* and *CACNA1C*, have provided two important insights into the nature of inter-individual risk for developing arrhythmias in response to drugs that block I_{Kr} , the major cause of drug-induced cardiac arrhythmias.¹⁹ First, cells with low G_{CaL} (and hence low G_{Kr} at baseline) exhibited the greatest prolongation of AP duration when exposed to I_{Kr} drug block (Figures 5 and 7). Second, cells with high G_{CaL} (and hence high G_{Kr} at baseline) showed greater propensity for development of EADs at moderate levels of I_{Kr} drug block (Figure 5). This observation also provides a plausible explanation for the well-described clinical observation that prolongation of the QT interval is an imperfect surrogate marker for proarrhythmic risk.⁴³ A corollary of the observation that a high G_{CaL} increases the susceptibility to EADs in response to drug block of I_{Kr} is that the co-

administration of an I_{CaL} blocker should reduce the risk of EADs (Supplementary material online, Figure S8). This is consistent with the observation that the administration of magnesium, which is a mild calcium channel blocker, is helpful in the acute management of patients with drug-induced *torsades de pointes*,¹⁹ and conversely that hypomagnesaemia, which would stimulate I_{CaL} , can exacerbate *torsades de pointes*.⁴⁴ It is also consistent with the observation that drugs that block I_{CaL} and I_{Kr} (e.g. verapamil) are not associated with drug-induced arrhythmias⁴⁵ and that verapamil prevented the development of *torsades de pointes* in rabbit hearts exposed to an I_{Kr} blocker.⁴⁶ However, given that calcium channel blockers are contra-indicated in some ventricular arrhythmias,⁴⁷ and the likelihood that patients who have drug-induced LQTS may have other underlying cardiac conditions,¹⁹ one should be cautious about prescribing calcium channel blockers. Conversely, it would be reasonable to consider using calcium channel blockers to treat patients with LQTS type 2 (i.e. patients with an isolated loss of I_{Kr} function) who continue to have cardiac events despite treatment with β -blockers.⁴⁸

4.1 Limitations

It is much more problematic to study gene expression in adult human samples, as opposed to laboratory animals, as there is necessarily a much longer time interval between death and obtaining tissue samples to process for mRNA expression. For this reason, in addition to analysing the full GTEx dataset, we also restricted our analysis to only those samples where the RIN value was >7 , which resulted in exclusion of \sim half of the samples. Similarly, it is difficult to obtain human heart samples for functional studies. For this reason, we used human iPSC studies. It should be noted, however, that some ion channels important for function in adult cardiac myocytes are not expressed at significant levels in immature cardiac myocytes⁴⁹ and so iPSC-CM may not be the ideal system for looking at ion channel modules other than *KCNH2* and *CACNA1C*. A further potential limitation of using iPSC-derived cardiac myocytes is the possibility that differences in maturity between wells and between cell lines could contribute to the differences we observed in our functional studies. This does not detract from the observation that wells that contained cells with higher levels of expression of *CACNA1C* and *KCNH2* showed less prolongation of FPD in response to I_{Kr} drug block, however, we acknowledge that the inter-individual differences we observed in immature cardiac myocytes may be larger than the differences one might see in adult myocytes.

4.2 Conclusion

Co-expression of *CACNA1C* and *KCNH2* ion channel genes in the heart enhances the robustness of cardiac electrical activity. Furthermore, we show that the subset of the population who have above average levels of calcium channel expression are most at risk from arrhythmias caused by drug block of the rapid delayed rectifier potassium current. We also suggest that the combination of technologies we have used should be applicable to identifying modules of co-expressed genes important in other fundamental physiological processes as well as investigating how the different modules of co-expressed genes interact to co-ordinate whole-of-cell function. Similar approaches should also be applicable to investigating the transcription factors and/or networks that control the co-expression of rhythmome gene modules.

Supplementary material

Supplementary material is available at *Cardiovascular Research* online.

Acknowledgements

This work was supported by the Victor Chang Cardiac Research Institute Innovation Centre, funded by the NSW Government. The Genotype-Tissue Expression (GTEx) Project was supported by the Common Fund of the Office of the Director of the National Institutes of Health, and by NCI, NHGRI, NHLBI, NIDA, NIMH, and NINDS. The data used for the analyses described in this article were obtained from the GTEx Portal on 16 February 2020.

Conflict of interest: none declared.

Funding

This work was supported by grants from the National Health and Medical Research Council (NHMRC) (App1116948 to J.I.V., App1074386 to J.I.V., and App1164518 to A.P.H.) and the National Institutes of Health (NIH) (R01LM012736 to J.A.G. and R01MH113005 to J.A.G.).

Data availability

Raw data for MEA recordings have been deposited at DRYAD database: doi: <https://doi.org/10.5061/dryad.612jm640k>. Details of the Nanostring codes and the raw data for mRNA expression levels have been deposited at National Center for Biotechnology Information Gene Expression Omnibus database under accession GSE146667. Code for the analysis of the public RNA Seq datasets is available at <https://github.com/sarbal/hERG-cal> and for the AP modelling is available at <https://git.victorchang.edu.au/projects/CC/repos/ordcoreg>.

References

- Zipes DP, Wellens HJ. Sudden cardiac death. *Circulation* 1998;**98**:2334–2351.
- Goyal V, Jassal DS, Dhalla NS. Pathophysiology and prevention of sudden cardiac death. *Can J Physiol Pharmacol* 2016;**94**:237–244.
- Moss AJ. MADIT-I and MADIT-II. *J Cardiovasc Electrophysiol* 2003;**14**(Suppl. 9): S96–S98.
- Behr E, Ensam B. New approaches to predicting the risk of sudden death. *Clin Med* 2016;**16**:283–283.
- Marder E, Goaillard J-M. Variability, compensation and homeostasis in neuron and network function. *Nat Rev Neurosci* 2006;**7**:563–574.
- Weiss JN, Karma A, MacLellan WR, Deng M, Rau CD, Rees CM, Wang J, Wisniewski N, Eskin E, Horvath S, Qu Z, Wang Y, Lusa AJ. “Good enough solutions” and the genetics of complex diseases. *Circ Res* 2012;**111**:493–504.
- Prinz AA, Billimoria CP, Marder E. Alternative to hand-tuning conductance-based models: construction and analysis of databases of model neurons. *J Neurophysiol* 2003;**90**:3998–4015.
- Britton OJ, Bueno-Orovio A, Van Ammel K, Lu HR, Towart R, Gallacher DJ, Rodriguez B. Experimentally calibrated population of models predicts and explains intersubject variability in cardiac cellular electrophysiology. *Proc Natl Acad Sci U S A* 2013;**110**:E2098–E2105.
- Sarkar AX, Sobie EA. Quantification of repolarization reserve to understand inter-patient variability in the response to pro-arrhythmic drugs: a computational analysis. *Heart Rhythm* 2011;**8**:1749–1755.
- Bányász T, Horváth B, Jian Z, Izu LT, Chen-Izu Y. Sequential dissection of multiple ionic currents in single cardiac myocytes under action potential-clamp. *J Mol Cell Cardiol* 2011;**50**:578–581.
- Rees CM, Yang J-H, Santolini M, Lusa AJ, Weiss JN, Karma A. The Ca²⁺ transient as a feedback sensor controlling cardiomyocyte ionic conductances in mouse populations. *eLife* 2018;**7**:4417.
- Huang CL-H. Murine electrophysiological models of cardiac arrhythmogenesis. *Physiol Rev* 2017;**97**:283–409.
- O’Leary T, Sutton AC, Marder E. Computational models in the age of large datasets. *Curr Opin Neurobiol* 2015;**32**:87–94.
- Schulz DJ, Goaillard J-M, Marder EE. Quantitative expression profiling of identified neurons reveals cell-specific constraints on highly variable levels of gene expression. *Proc Natl Acad Sci U S A* 2007;**104**:13187–13191.
- Eisen MB, Spellman PT, Brown PO, Botstein D. Cluster analysis and display of genome-wide expression patterns. *Proc Natl Acad Sci U S A* 1998;**95**:14863–14868.

- Gaiteri C, Ding Y, French B, Tseng GC, Sibille E. Beyond modules and hubs: the potential of gene coexpression networks for investigating molecular mechanisms of complex brain disorders. *Genes Brain Behav* 2014;**13**:13–24.
- Lee HK, Hsu AK, Sajdak J, Qin J, Pavlidis P. Coexpression analysis of human genes across many microarray data sets. *Genome Res* 2004;**14**:1085–1094.
- Ballouz S, Verleyen W, Gillis J. Guidance for RNA-seq co-expression network construction and analysis: safety in numbers. *Bioinformatics* 2015;**31**:2123–2130.
- Roden DM. Drug-induced prolongation of the QT interval. *N Engl J Med* 2004;**350**:1013–1022.
- Zoubarev A, Hamer KM, Keshav KD, McCarthy EL, Santos JRC, Van Rossum T, McDonald C, Hall A, Wan X, Lim R, Gillis J, Pavlidis P. Gemma: a resource for the re-use, sharing and meta-analysis of expression profiling data. *Bioinformatics* 2012;**28**:2272–2273.
- The GTEx Consortium. The Genotype-Tissue Expression (GTEx) project. *Nat Genet* 2013;**45**:580–585.
- Burridge PV, Diecke S, Matsa E, Sharma A, Wu H, Wu JC. Modeling cardiovascular diseases with patient-specific human pluripotent stem cell-derived cardiomyocytes. *Methods Mol Biol* 2016;**1353**:119–130.
- Mills RJ, Titmarsh DM, Koenig X, Parker BL, Ryall JG, Quaife-Ryan GA, Voges HK, Hodson MP, Ferguson C, Drowley L, Plowright AT, Needham EJ, Wang QD, Gregorevic P, Xin M, Thomas WG, Parton RG, Nielsen LK, Launikonis BS, James DE, Elliott DA, Porrello ER, Hudson JE. Functional screening in human cardiac organoids reveals a metabolic mechanism for cardiomyocyte cell cycle arrest. *Proc Natl Acad Sci U S A* 2017;**114**:E8372–E8381.
- Perry MD, Ng CA, Mangala MM, Ng TYM, Hines AD, Liang W, Xu MJO, Hill AP, Vandenberg JL. Pharmacological activation of I_{Kr} in models of long QT Type 2 risks overcorrection of repolarization. *Cardiovasc Res* 2020;**116**:1434–1445.
- O’Hara T, Virág L, Varró A, Rudy Y. Simulation of the un diseased human cardiac ventricular action potential: model formulation and experimental validation. *PLoS Comput Biol* 2011;**7**:e1002061.
- Mann SA, Intiaz M, Winbo A, Rydberg A, Perry MD, Couderc JP, Polonsky B, McNitt S, Zareba W, Hill AP, Vandenberg JL. Convergence of models of human ventricular myocyte electrophysiology after global optimization to recapitulate clinical long QT phenotypes. *J Mol Cell Cardiol* 2016;**100**:25–34.
- Krogh-Madsen T, Jacobson AF, Ortega FA, Christini DJ. Global optimization of ventricular myocyte model to multi-variable objective improves predictions of drug-induced torsades de pointes. *Front Physiol* 2017;**8**:e1000173.
- Heitmann S, Aburn MJ, Breakspear M. The brain dynamics toolbox for Matlab. *Neurocomputing* 2018;**315**:82–88.
- Sadrieh A, Mann SA, Subbiah RN, Domanski L, Taylor JA, Vandenberg JL, Hill AP. Quantifying the origins of population variability in cardiac electrical activity through sensitivity analysis of the electrocardiogram. *J Physiol* 2013;**591**:4207–4222.
- Bush WS, Crawford DC, Alexander C, George AL Jr, Roden DM, Ritchie MD. Genetic variation in the rhythmome: ethnic variation and haplotype structure in candidate genes for arrhythmias. *Pharmacogenomics* 2009;**10**:1043–1053.
- Sah R, Ramirez RJ, Backx PH. Modulation of Ca²⁺ release in cardiac myocytes by changes in repolarization rate. *Circ Res* 2002;**90**:165–173.
- Shimizu W, Moss AJ, Wilde AAM, Towbin JA, Ackerman MJ, January CT, Tester DJ, Zareba W, Robinson JL, Qi M, Vincent GM, Kaufman ES, Hofman N, Noda T, Kamakura S, Miyamoto Y, Shah S, Amin V, Goldenberg I, Andrews ML, McNitt S. Genotype-phenotype aspects of type 2 long QT syndrome. *J Am Coll Cardiol* 2009;**54**:2052–2062.
- Harris K, Aylott M, Cui Y, Louttit JB, McMahon NC, Sridhar A. Comparison of electrophysiological data from human-induced pluripotent stem cell-derived cardiomyocytes to functional preclinical safety assays. *Toxicol Sci* 2013;**134**:412–426.
- Pedersen HS, Elming H, Seibæk M, Burchardt H, Brendorp B, Torp-Pedersen C, Køber L. Risk factors and predictors of Torsade de Pointes ventricular tachycardia in patients with left ventricular systolic dysfunction receiving dofetilide. *Am J Cardiol* 2007;**100**:876–880.
- Sayed N, Liu C, Wu JC. Translation of human-induced pluripotent stem cells. *J Am Coll Cardiol* 2016;**67**:2161–2176.
- Eichel CA, Ríos-Pérez EB, Liu F, Jameson MB, Jones DK, Knickebine JJ, Robertson GA. A microtranslatome coordinately regulates sodium and potassium currents in the human heart. *Elife* 2019;**8**:e52654.
- Min B, Zheng M. Correlated network of networks enhances robustness against catastrophic failures. *PLoS One* 2018;**13**:e0195539.
- Lundby A, Rossin EJ, Steffensen AB, Acha MR, Newton-Cheh C, Pfeufer A, Lynch SN, QT Interval International GWAS Consortium (QT-IGC), Olesen SP, Brunak S, Ellinor PT, Jukema JW, Trompet S, Ford I, Macfarlane PW, Krijthe BP, Hofman A, Uitterlinden AG, Stricker BH, Nathoe HM, Spiering W, Daly MJ, Asselbergs FW, van der Harst P, Milan DJ, de Bakker PI, Lage K, Olsen JV. Annotation of loci from genome-wide association studies using tissue-specific quantitative interaction proteomics. *Nat Methods* 2014;**11**:868–874.
- Nerbonne JM, Kass RS. Molecular physiology of cardiac repolarization. *Physiol Rev* 2005;**85**:1205–1253.
- Ponce-Balbuena D, Guerrero-Serna G, Valdivia CR, Caballero R, Diez-Guerra FJ, Jiménez-Vázquez EN, Ramírez RJ, Monteiro da Rocha A, Herron TJ, Campbell KF,

- Willis BC, Alvarado FJ, Zarzoso M, Kaur K, Pérez-Hernández M, Matamoros M, Valdivia HH, Delpón E, Jalife J. Cardiac Kir2.1 and Nav1.5 channels traffic together to the sarcolemma to control excitability. *Circ Res* 2018;**122**:1501–1516.
41. Milstein ML, Musa H, Balbuena DP, Anumonwo JM, Auerbach DS, Fursan PB, Hou L, Hu B, Schumacher SM, Vaidyanathan R, Martens JR, Jalife J. Dynamic reciprocity of sodium and potassium channel expression in a macromolecular complex controls cardiac excitability and arrhythmia. *Proc Natl Acad Sci U S A* 2012;**109**:E2134–E2143.
42. Deschênes I, Armoundas AA, Jones SP, Tomaselli GF. Post-transcriptional gene silencing of KChIP2 and Navbeta1 in neonatal rat cardiac myocytes reveals a functional association between Na and Ito currents. *J Mol Cell Cardiol* 2008;**45**:336–346.
43. Sager PT, Gintant G, Turner JR, Pettit S, Stockbridge N. Rechanneling the cardiac proarrhythmia safety paradigm: a meeting report from the Cardiac Safety Research Consortium. *Am Heart J* 2014;**167**:292–300.
44. Roden DM, Iansmith DH. Effects of low potassium or magnesium concentrations on isolated cardiac tissue. *Am J Med* 1987;**82**:18–23.
45. Redfern WS, Carlsson L, Davis AS, Lynch WG, MacKenzie I, Palethorpe S, Siegl PK, Strang I, Sullivan AT, Wallis R, Camm AJ, Hammond TG. Relationships between pre-clinical cardiac electrophysiology, clinical QT interval prolongation and torsade de pointes for a broad range of drugs: evidence for a provisional safety margin in drug development. *Cardiovasc Res* 2003;**58**:32–45.
46. Farkas AS, Makra P, Csik N, Orosz S, Shattock MJ, Fülöp F, Forster T, Csanády M, Papp JG, Varró A, Farkas A. The role of the Na⁺/Ca²⁺ exchanger, INa and ICaL in the genesis of dofetilide-induced torsades de pointes in isolated, AV-blocked rabbit hearts. *Br J Pharmacol* 2009;**156**:920–932.
47. Akhtar M, Tchou P, Jazayeri M. Use of calcium channel entry blockers in the treatment of cardiac arrhythmias. *Circ* 1989;**80**:IV31–IV39.
48. Komiya N, Tanaka K, Doi Y, Fukae S, Nakao K, Isomoto S, Seto S, Yano K. A patient with LQTS in whom verapamil administration and permanent pacemaker implantation were useful for preventing torsade de pointes. *Pacing Clin Electrophysiol* 2004;**27**:123–124.
49. Robertson C, Tran DD, George SC. Concise review: maturation phases of human pluripotent stem cell-derived cardiomyocytes. *Stem Cells* 2013;**31**:829–837.

Translational perspective

Here, we show, using meta-analysis of multiple independent human gene expression datasets, that there is co-expression of *KCNH2*–*CACNA1C* in human heart tissue which was then confirmed in human cardiac myocytes derived from induced pluripotent stem cells. Both *in silico* and functional studies show that the co-expression of *CACNA1C* and *KCNH2* increases the robustness of cardiac electrical signalling. Our data also suggest that those patients who express higher levels of *CACNA1C* and *KCNH2* are likely to be more susceptible to arrhythmias when exposed to drugs that block I_{Kr} , the major cause of drug-induced cardiac arrhythmias.