

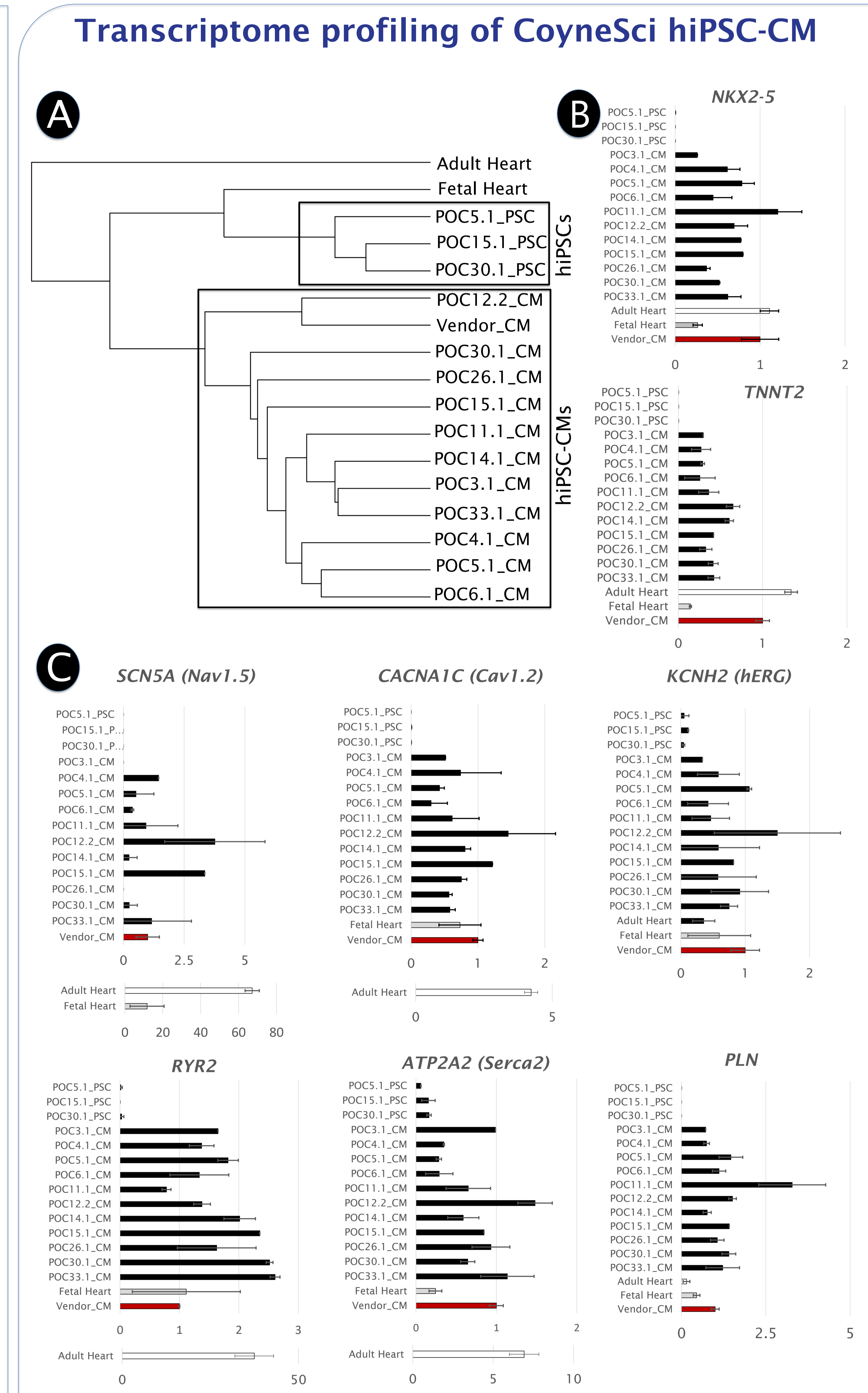
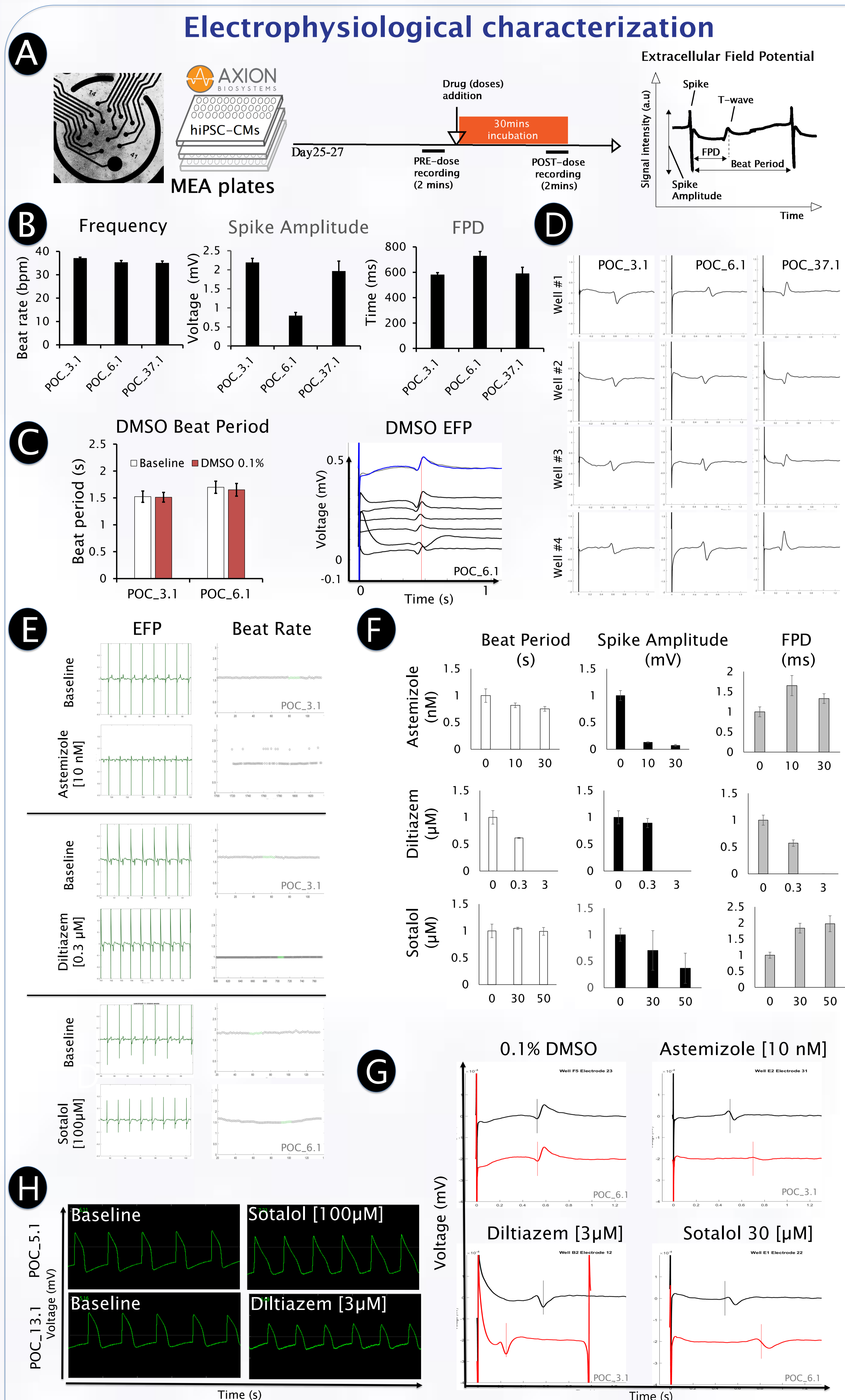


Inter-Individual Differences in Electrophysiology, Pharmacology and Drug Responses in a Cohort of hiPSC-derived Cardiomyocytes: Endorsing Clinical Trials-in-a-Dish

Bernard Fermini, Denise Sullivan, Jerome Chal, Eli Fine, Alexander Ulan Hernandez, Alexia King, Jessica Johnson, Shawn Coyne and Kevin P. Coyne
Coyne Scientific, Atlanta, GA, USA

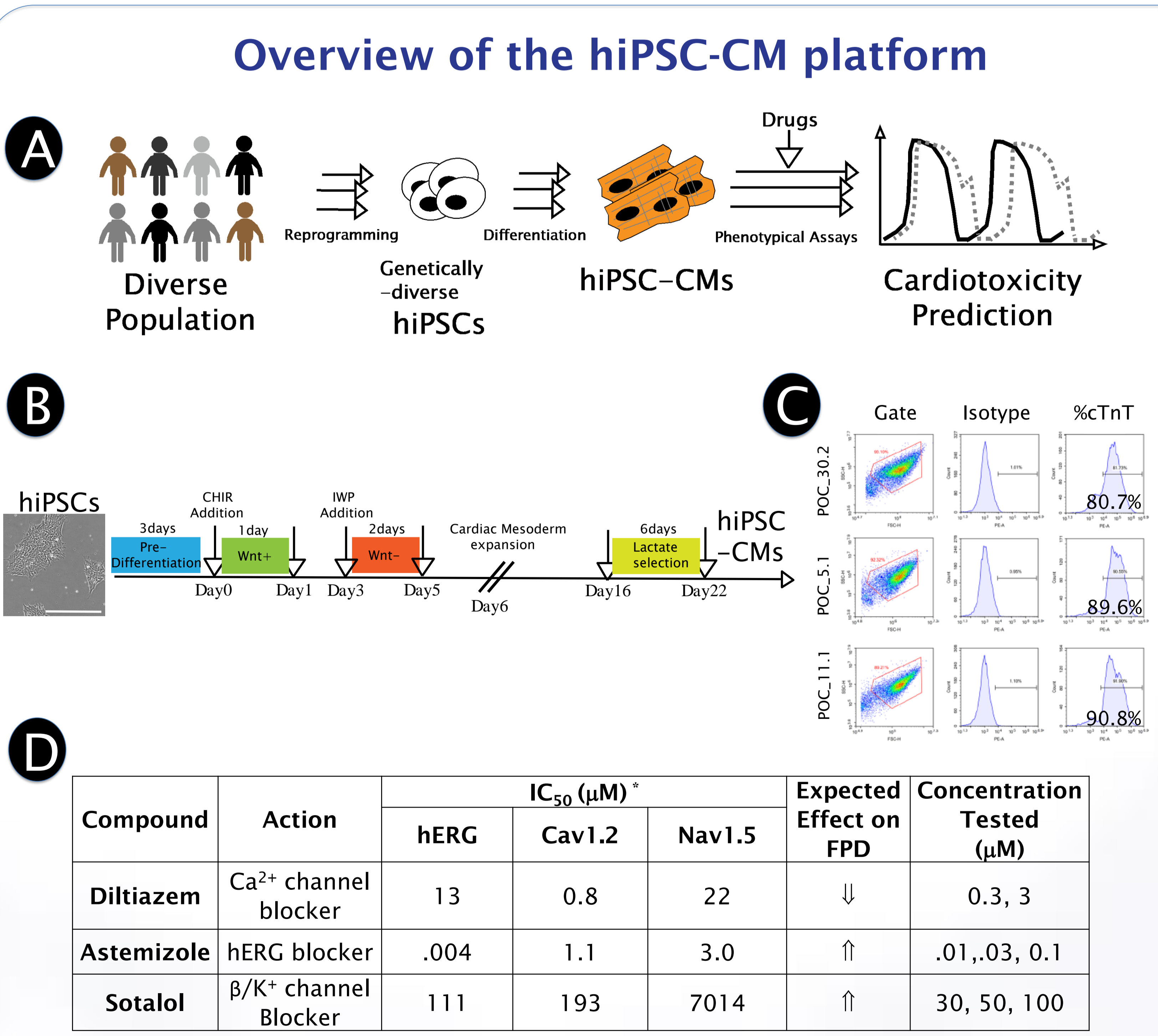


The concept of “Clinical trials-in-a-dish” or *in vitro* clinical trials, has received significant attention lately because of emerging technologies that allow testing novel compounds on patient cells before moving into actual clinical trials. Moreover, recent studies have shown that individual serious adverse events (SAEs) susceptibility in a population of volunteers with unknown genetic background can be recapitulated in human inducible pluripotent stem cell cardiomyocytes (hiPSC-CMs) derived from these same individuals, providing a proof of concept for *in vitro* preclinical trials. Hence, leveraging genetic diversity in preclinical testing may hold the key to reducing clinical attrition, based on an increased ability to predict SAEs in a population. In this study, we examined the electrophysiological and pharmacological properties of hiPSC-CMs derived from healthy donors. We first established the full transcriptome profiling of our cell lines, and found inter-line differences in the expression levels of important cardiac markers, and key ion channels. We then defined individual electrophysiological and pharmacological properties of a number of lines by measuring various parameters extracted from extracellular field potential (EFP) and Local Extracellular Action Potential (LEAP) recordings from spontaneously beating cells. Finally, we studied the effects of 3 different CiPA drugs (from all 3 risk categories for arrhythmia liability), on our hiPSC-CMs to establish a pharmacological profile.



(A) Whole transcriptome analysis of hiPSCs (PSC) and hiPSC-CMs in comparison to commercially available hiPSC-CMs (Vendor_CM) and human heart tissues. Data were generated by TempO-Seq (BioSpyder) whole transcriptome assay. Dendrogram was generated by maximum pairwise difference clustering. (B) Comparison of expression of cardiac markers and (C) various ion channels and regulators between hiPSCs and hiPSC-CMs. Values are normalized to Vendor CM sample.

Conclusions
Our data demonstrate that hiPSC-CMs derived from healthy donors exhibit electrophysiological and pharmacological profiles that are unique to each donor, resulting in a distribution of drug effects and potencies. Our results show line-specific baseline properties supporting validation and implementation of testing for safety and toxicity of drugs across a cohort of donors, rather than a single cell line, endorsing the concept of Clinical Trials-in-a-Dish.



*From: Kramer, J.; Obejero-Paz, C. A.; Myatt, G. et al., MICE Models: Superior to the HERG Model in Predicting Torsades de Pointes. *Sci. Rep.* 2013, 3. FPD: Field potential duration.

(A) Cohort of hiPSCs was generated from a genetically-diverse pool of donors, differentiated into hiPSC-CMs, and used to define their electrophysiological and pharmacological profile. (B) Directed differentiation protocol for hiPSC-CMs production, adapted from the GiWi protocol of Palecek and colleagues. (C) Representative flow cytometric analysis of cTnT for 3 lines of hiPSC-CMs. (D) Table of compound descriptions and concentrations used to dose hiPSC-CMs.