

# Robust and Efficient Generation of Functional Human Pluripotent Stem Cell-Derived Atrial Cardiomyocytes

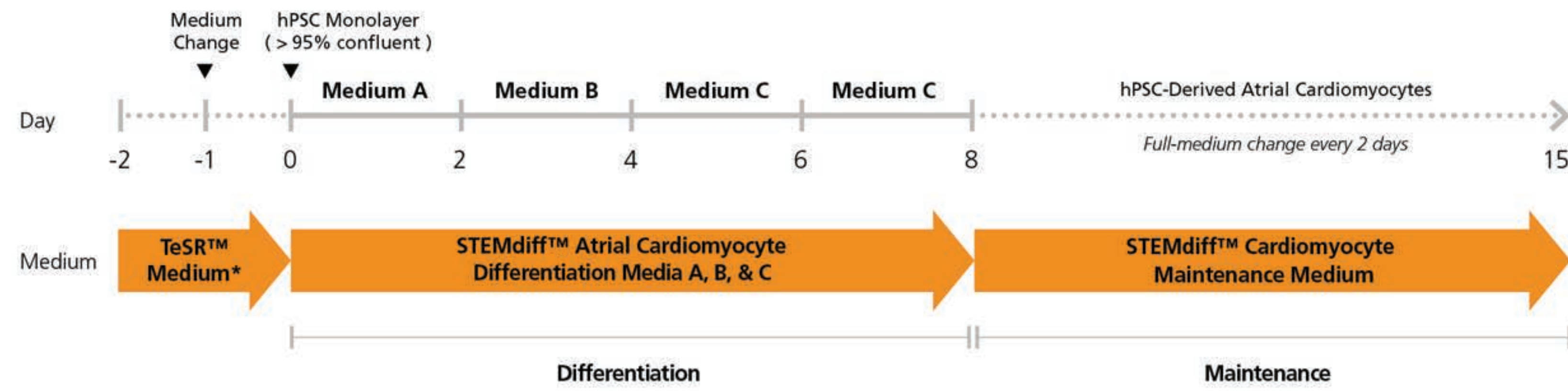
Jessica Legree<sup>1</sup>, Trisha Ang<sup>1</sup>, Vivian Nguyen<sup>1</sup>, Hattie Luo<sup>1</sup>, Steve S. Szilvassy<sup>1</sup>, Allen C. Eaves<sup>1,2</sup>, Sharon A. Louis<sup>1</sup>, and Vincenzo Macri<sup>1</sup>

<sup>1</sup>STEMCELL Technologies Inc., Vancouver BC, Canada; <sup>2</sup>Terry Fox Laboratory, BC Cancer, Vancouver BC, Canada

## INTRODUCTION

The ability to generate and maintain high-quality human pluripotent stem cell (hPSC)-derived atrial cardiomyocytes is necessary to model atrial cardiomyocyte disease and for drug discovery and toxicology. Protocols and reagents to make hPSC-derived atrial cardiomyocytes are variable and not standardized. We have developed the STEMdiff™ Atrial Cardiomyocyte Differentiation Kit, a defined, serum-free cell culture medium to differentiate hPSCs to atrial cardiomyocytes. This kit provides a standardized differentiation workflow to generate functional hPSC-derived atrial cardiomyocytes that are ready for use in downstream applications.

## METHODS

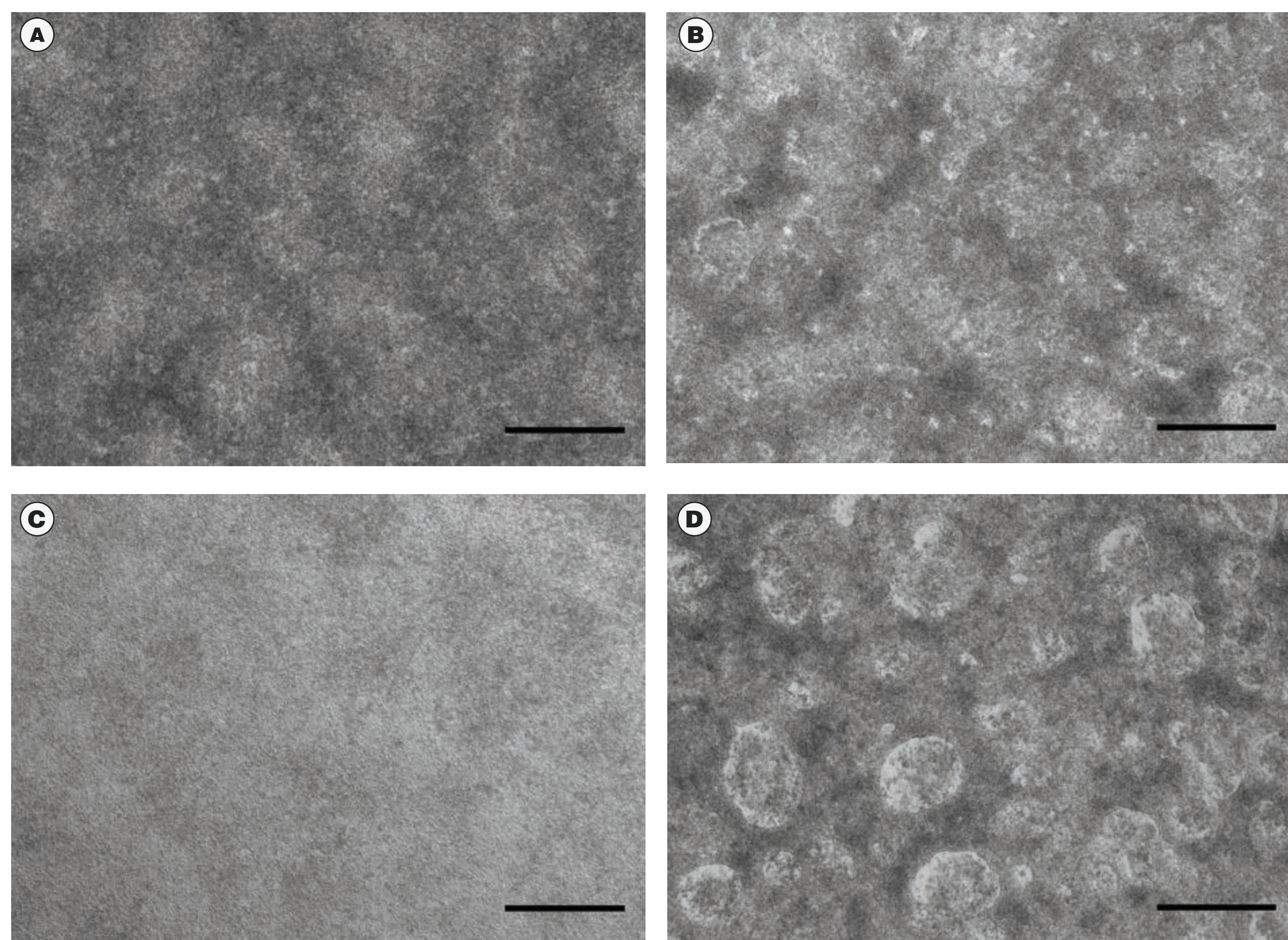


\*mTeSR™1, mTeSR™ Plus, TeSR™-AOF, or TeSR™-E8™

**FIGURE 1. STEMdiff™ Atrial Cardiomyocyte Differentiation Kit Workflow**

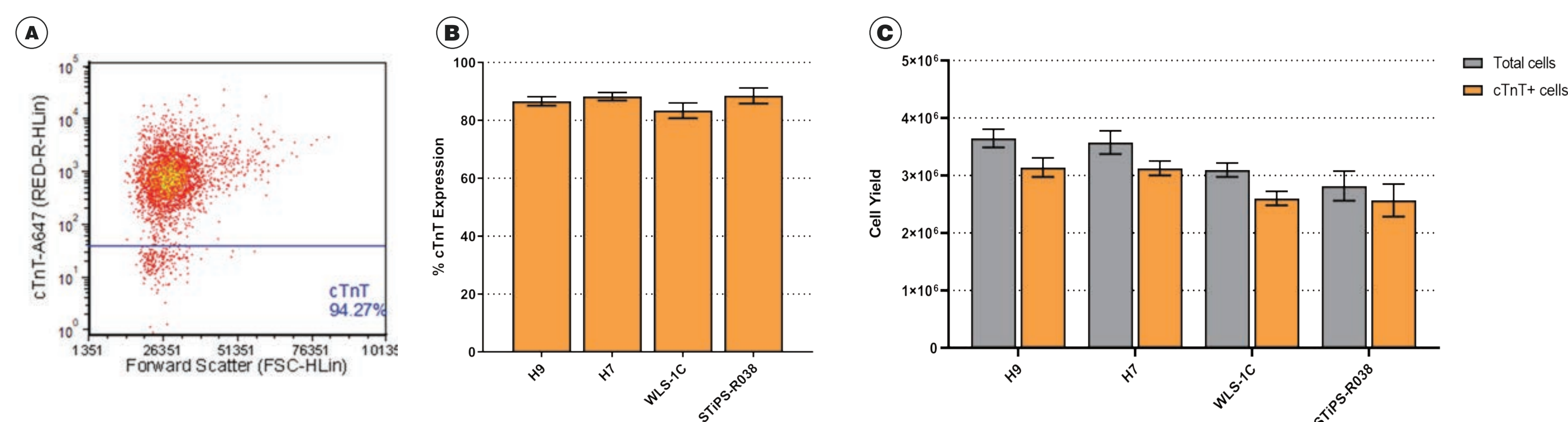
hPSCs were seeded as single cells ( $3.5 - 8 \times 10^5$  cells/well) onto Matrigel®-coated 12-well plates in TeSR™ medium containing Y-27632 and maintained for two days. On day 0, differentiation of hPSCs to atrial cardiomyocytes was initiated with a full-medium change using Differentiation Medium A. On days 2 and 4, the medium was replaced with Differentiation Medium B and C, respectively. On day 8, Differentiation Medium C was replaced with STEMdiff™ Cardiomyocyte Maintenance Medium. A full-medium change was completed every other day until day 15. On day 15, a confluent, beating monolayer was observed, and the hPSC-derived atrial cardiomyocytes were harvested for characterization using flow cytometry, qPCR, RNAseq, and electrophysiology.

## RESULTS



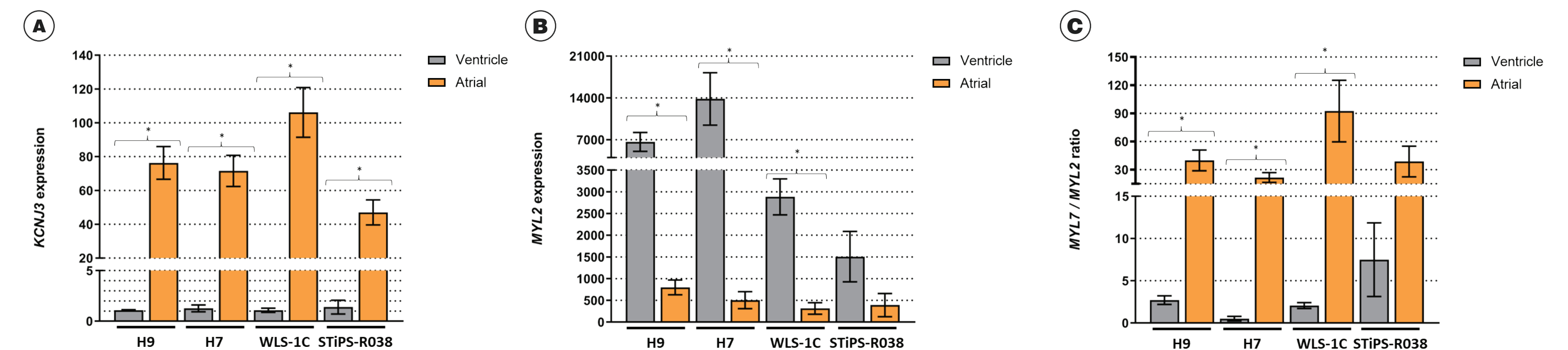
**FIGURE 2. STEMdiff™ Atrial Cardiomyocyte Differentiation Kit Produced Robust, Uniform Monolayers of hPSC-derived Atrial Cardiomyocytes Across hPSC Lines**

hPSC-derived atrial cardiomyocytes generated by the STEMdiff™ Atrial Cardiomyocyte Differentiation Kit exhibited confluent, beating monolayers on day 15 of differentiation. Robust, uniform beating monolayers were observed in four hPSC lines, including 2 embryonic stem (ES) cell lines (A) H9 and (B) H7 and 2 induced pluripotent stem (iPS) cell lines (C) WLS-1C and (D) STiPS-R038. Scale bar = 500 μm.



**FIGURE 3. STEMdiff™ Atrial Cardiomyocyte Differentiation Kit Generated a High Percentage and Yield of cTnT-Positive Cells Across hPSC Lines**

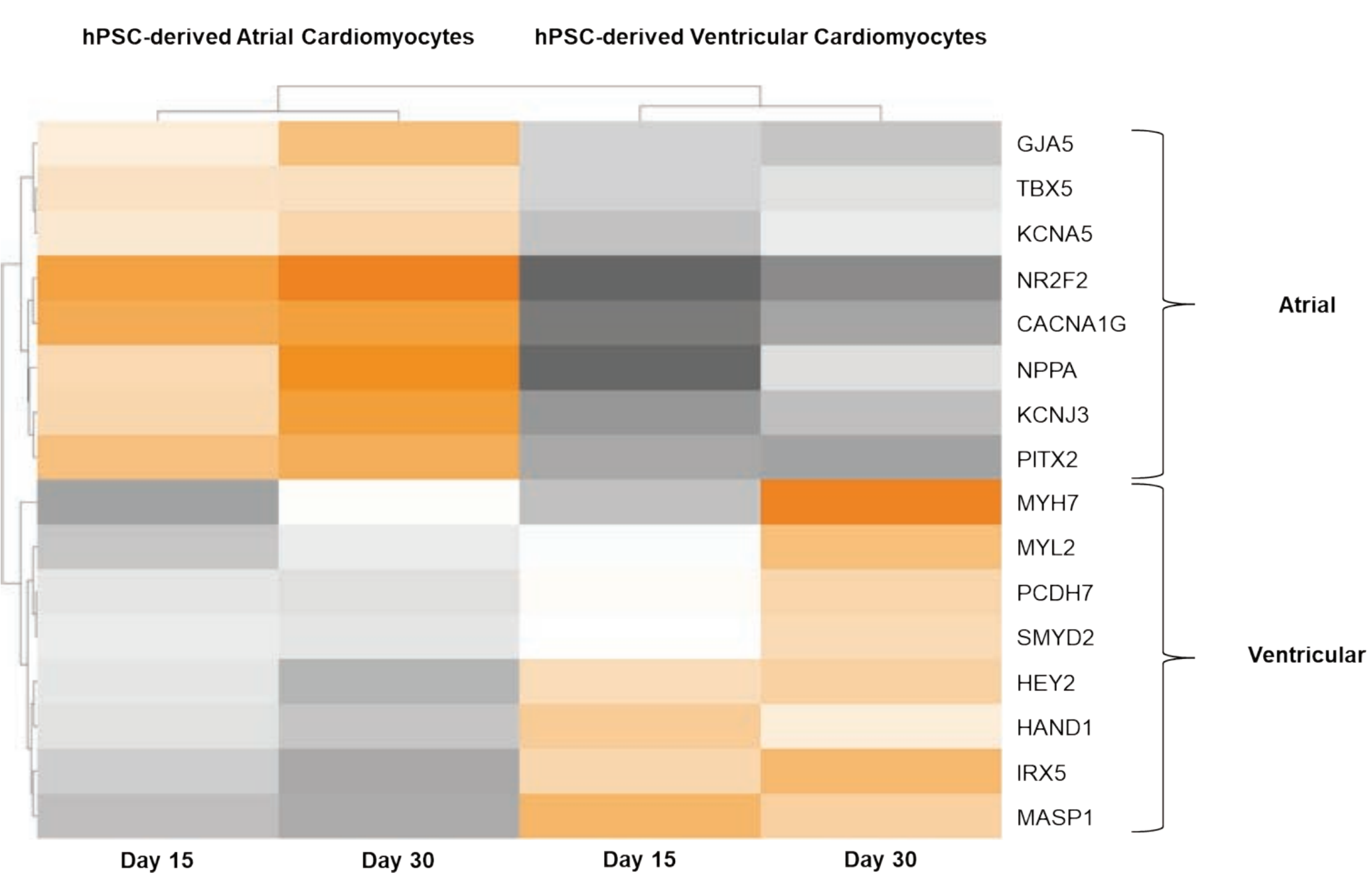
To demonstrate the efficiency of the STEMdiff™ Atrial Cardiomyocyte Differentiation Kit, four hPSC lines were differentiated to atrial cardiomyocytes and were harvested for flow cytometry. Differentiation was performed in a 12-well plate format. (A) Flow cytometry histogram of cTnT expression for cardiomyocytes generated with the STiPS-R038 hiPSC line. Bar graphs showing (B) > 80% cTnT expression and (C) > 2 million cardiomyocytes were produced from four hPSC lines. Data are shown as mean ± SEM, n = 18.



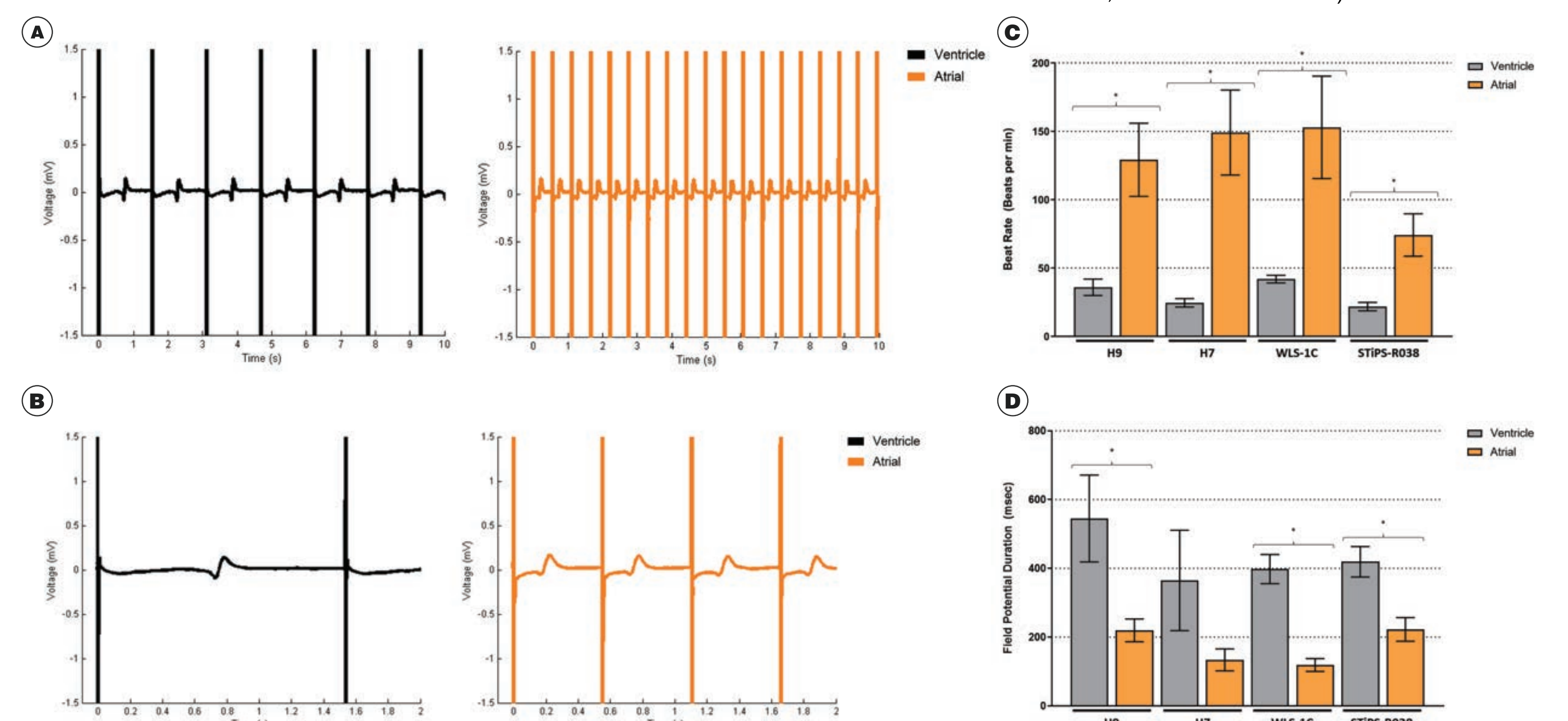
**FIGURE 4. Quantitative PCR Revealed Upregulation of *KCNJ3* and Downregulation of *MYL2* in hPSC-derived Atrial Cardiomyocytes**

Four hPSC lines were differentiated to either atrial cardiomyocytes using the STEMdiff™ Atrial Cardiomyocyte Differentiation Kit or ventricular cardiomyocytes using the STEMdiff™ Ventricular Cardiomyocyte Differentiation Kit. Cardiomyocyte subtypes were assessed by qPCR for expression of (A) *KCNJ3*, an atrial-specific potassium ion channel gene, and (B) *MYL2*, a ventricle-specific regulatory light chain of myosin gene. (C) The ratio of *MYL7/MYL2* expression can be used as an indicator for an atrial-like gene expression. Across four hPSC lines, atrial cardiomyocytes showed significantly increased *KCNJ3* expression (> 30-fold, n = 40), significantly decreased *MYL2* expression (< 0.3-fold, n = 63), and significantly increased *MYL7/MYL2* ratio (> 2-fold, n = 63) compared to ventricular cardiomyocytes. Data are shown as mean ± SEM. Asterisk indicates significance of p < 0.05.

**FIGURE 5. RNAseq Timepoint Analysis Revealed Upregulation of Atrial-Specific Genes and Downregulation of Ventricle-Specific Genes in hPSC-Derived Atrial Cardiomyocytes**



Four hPSC lines were differentiated to either atrial cardiomyocytes using the STEMdiff™ Atrial Cardiomyocyte Differentiation Kit or ventricular cardiomyocytes using the STEMdiff™ Ventricular Cardiomyocyte Differentiation Kit. Cardiomyocyte subtypes were harvested on day 15 and day 30 for RNAseq analyses. hPSC-derived atrial cardiomyocytes showed upregulation of atrial-specific genes (*GJA5*, *TBX5*, *KCNA5*, *NR2F2*, *CACNA1G*, *NPPA*, *KCNJ3*, *PITX2*) and downregulation of ventricle-specific genes (*MYH7*, *MYL2*, *PCDH7*, *SMYD2*, *HEY2*, *HAND1*, *IRX5*, *MASP1*) on day 15 and day 30 compared to hPSC-derived ventricular cardiomyocytes. RNAseq data was generated from 3 hPSC lines (H9, WLS-1C, and STiPS-M001).



**FIGURE 6. Microelectrode Array of hPSC-Derived Atrial Cardiomyocytes Showed Increased Beat Rate and Shorter Field Potential Duration Compared to hPSC-Derived Ventricular Cardiomyocytes**

Microelectrode Array (Maestro MEA, Axion BioSystems) recordings of hPSC-derived atrial cardiomyocytes (orange) and hPSC-derived ventricular cardiomyocytes (black). hPSC-derived atrial cardiomyocytes showed (A) increased beat rate and (B) decreased field potential duration (FPD) compared to hPSC-derived ventricular cardiomyocytes. (C) Beat rate recordings from 4 hPSC lines demonstrated that hPSC-derived atrial cardiomyocytes consistently beat faster compared to hPSC-derived ventricular cardiomyocytes. Across 4 cell lines, the average atrial beat rate was  $123 \pm 14$  beats per minute (mean ± SEM; n = 21), and the average ventricular beat rate was  $32 \pm 3$  beats per minute (mean ± SEM; n = 16). (D) Across 4 cell lines, FPD values were consistently shorter in hPSC-derived atrial cardiomyocytes compared to ventricular cardiomyocytes. Average atrial FPD was  $181 \pm 18$  msec (mean ± SEM; n = 21), and average ventricular FPD was  $443 \pm 49$  msec (mean ± SEM; n = 16). Asterisk indicates a significance of p < 0.05.

## Summary

- STEMdiff™ Atrial Cardiomyocyte Differentiation Kit is a standardized, defined, and serum-free differentiation workflow for simple and efficient generation of functional hPSC-derived atrial cardiomyocytes.
- hPSC-derived atrial cardiomyocytes generated using the STEMdiff™ Atrial Cardiomyocyte Differentiation Kit produced uniform beating monolayers and generated a high percentage and yield of cTnT-positive cells across hPSC lines.
- hPSC-derived atrial cardiomyocytes demonstrated upregulation of atrial-specific genes (*KCNJ3*, *GJA5*, *NPPA*, *PITX2*) and downregulation of ventricle-specific genes (*MYL2*, *MYH7*, *HEY2*, *HAND1*).
- Electrophysiology profile of hPSC-derived atrial cardiomyocytes showed increased beat rate and shorter field potential duration (FPD) compared to hPSC-derived ventricular cardiomyocytes.



Jessica Legree,  
Technical Scientist  
Research & Development  
STEMCELL Technologies  
jessica.legree@stemcell.com