

Efficient Production and Processing of Cardiomyocytes from Human Pluripotent Stem Cells Using STEMdiff™ Cardiomyocyte Products

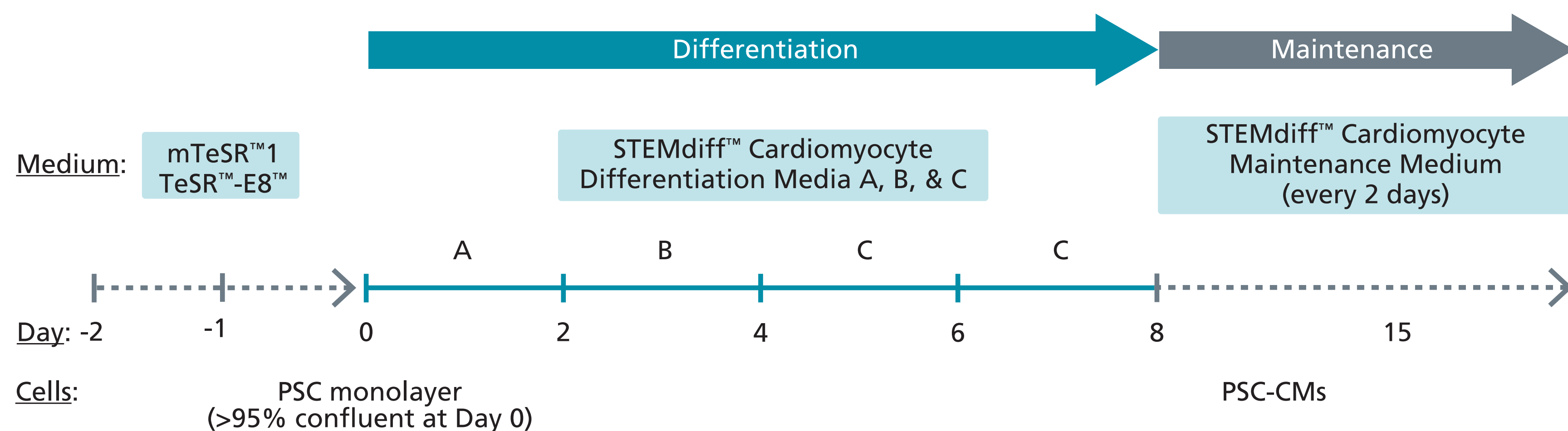
Vincenzo Macri¹, Jessica Norberg¹, Colleen Umali¹, Arthur V. Sampaio¹, Terry E. Thomas¹, Allen C. Eaves^{1,2}, Stephen J. Szilvassy¹, and Sharon Louis¹

¹STEMCELL Technologies Inc., Vancouver BC, Canada ²Terry Fox Laboratory, BC Cancer Agency, Vancouver BC, Canada

Introduction

Human pluripotent stem cell-derived cardiomyocytes (hPSC-CMs) are used for disease modelling, drug discovery, and toxicology screening. However, the efficiency of producing hPSC-CMs is variable and, once established in culture, processing of hPSC-CMs for downstream assays is cumbersome. To overcome these limitations, we have developed a complete workflow comprising several novel products: 1) STEMdiff™ Cardiomyocyte Differentiation Kit, 2) STEMdiff™ Cardiomyocyte Maintenance Kit, 3) STEMdiff™ Cardiomyocyte Dissociation Kit, 4) STEMdiff™ Cardiomyocyte Freezing Medium, and 5) STEMdiff™ Cardiomyocyte Support Medium.

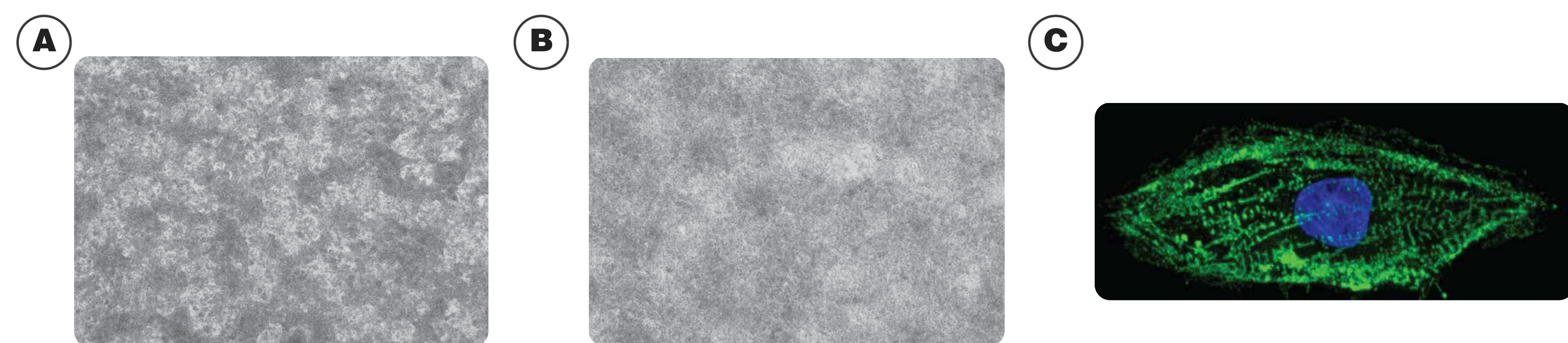
Methods



hPSCs were seeded as single cells (~3 x 10⁵/well) with Y-27632 ROCK inhibitor in TeSR™-E8™ or mTeSR™1 medium on Matrigel®-coated 12-well plates and maintained with daily medium changes until a confluent monolayer (>95%) was achieved. Day 0, the cells were treated with differentiation medium A. Day 2, medium A was removed and replaced with differentiation medium B. Day 4, medium B was removed and replaced with differentiation medium C. Day 6, medium C was removed and replaced with fresh differentiation medium C. Day 8, medium C was removed and replaced with maintenance medium. Maintenance medium was removed and replaced with fresh maintenance medium every 2 days until day 15. At Day 15, cardiomyocytes can be maintained up to a month or longer using the STEMdiff™ Cardiomyocyte Maintenance Kit or harvested with the STEMdiff™ Cardiomyocyte Dissociation Kit to replat for downstream applications (e.g. flow cytometry, immunocytochemistry, cryopreservation, electrophysiology, etc).

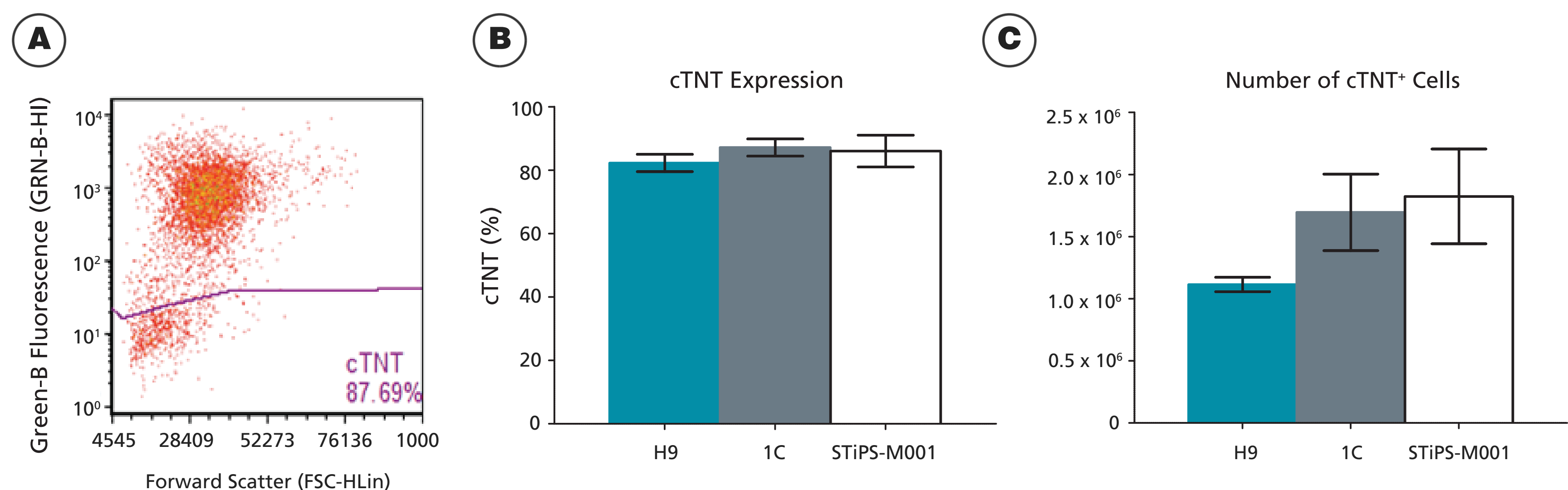
Results

FIGURE 1. hPSC-CMs generated with STEMdiff™ Cardiomyocyte Differentiation and Maintenance Kits



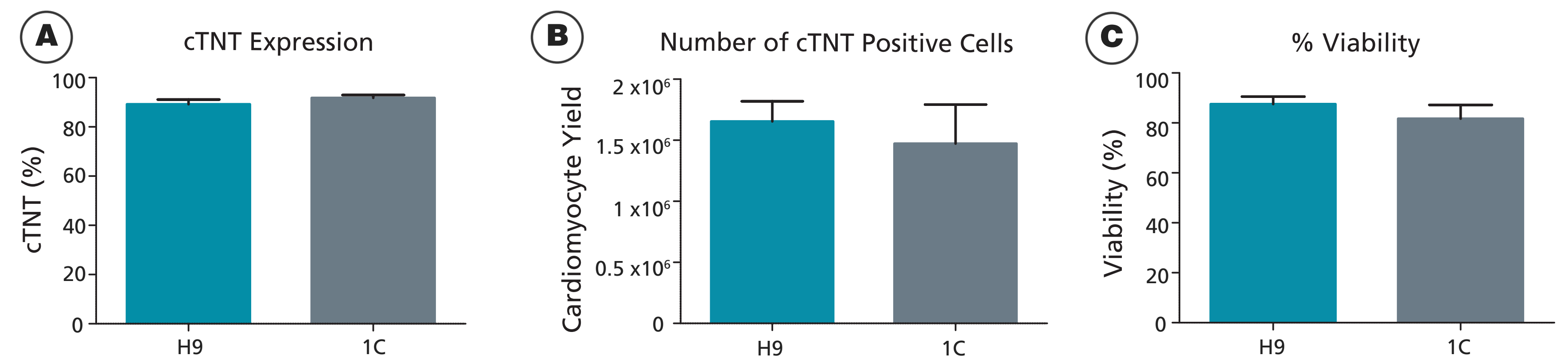
(A, B) Bright field images of hESC-CM and hiPSC-CM monolayers, respectively. (C) Confocal image of a single hiPSC-CM stained with cTNT (green) and Hoechst (blue).

FIGURE 2. STEMdiff™ Cardiomyocyte Differentiation and Maintenance Kits produce high percentage and yield of cTNT-positive cells from multiple human iPS and ES lines



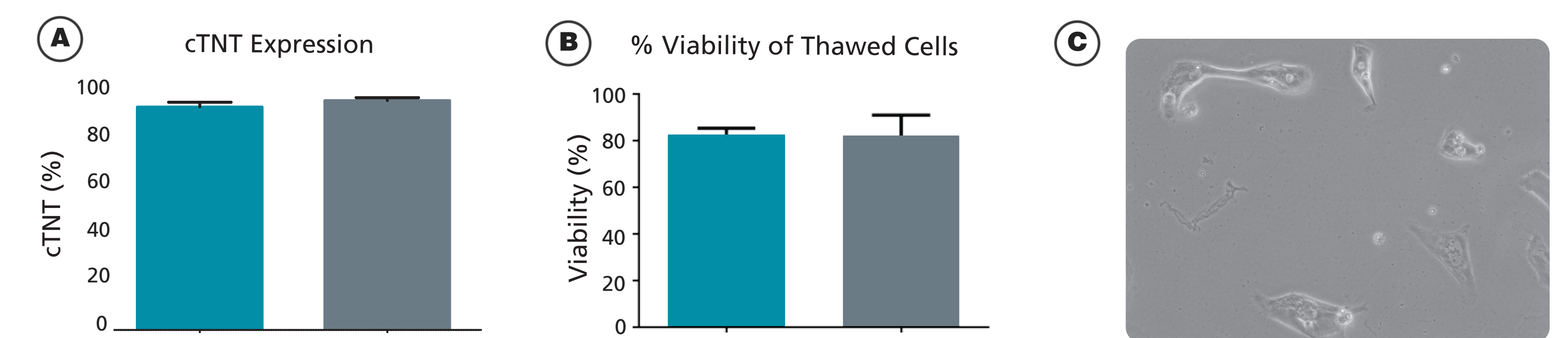
(A) Flow cytometry histogram of cTnT data for hPSC-CMs generated with the STiPS-M001 iPS line. (B, C) Bar graphs showing >80% cTNT-positive cells and >1 million cardiomyocytes are produced from hESC (H9) and hiPSC (1C, STiPS-M001) lines.

FIGURE 3. STEMdiff™ Cardiomyocyte Dissociation Kit harvests high yield of viable cardiomyocytes



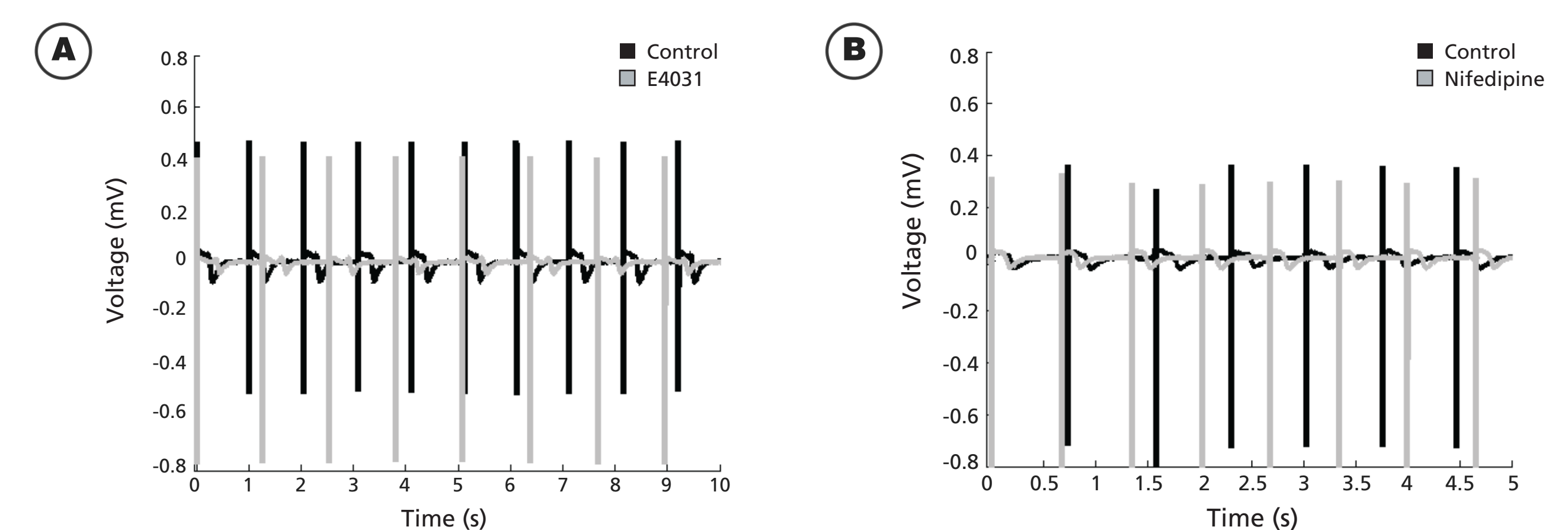
(A, B & C) hPSC-CMs have high purity (>85% cTnT-positive cells), yield (>1.5 million cardiomyocytes) and viability (>80% live cells) using the STEMdiff™ Cardiomyocyte Dissociation Kit. These values were determined from hPSC-CMs dissociated from a single well of a 12-well plate seeded initially with ES (H9) and iPS (1C) cells (n = 3).

FIGURE 4. STEMdiff™ Cardiomyocyte Freezing and Support Media used together support the cryopreservation, thawing and the replating of viable cardiomyocytes



(A, B) Cryopreserved hPSC-CMs have high purity (>85% cTnT-positive cells) and viability (>75% live cells) using the STEMdiff™ Cardiomyocyte Freezing Medium and after thawing with STEMdiff™ Cardiomyocyte Support Medium. These values were determined using 0.5 million cryopreserved hPSC-CMs at 1 and 3 months using ES (H9) and iPS (1C) cells (n = 2). (C) Bright field image of replated single hPSC-CMs at day 7 that were cryopreserved with STEMdiff™ Cardiomyocyte Freezing Medium and after thawing with STEMdiff™ Cardiomyocyte Support Medium. The cryopreserved hPSC-CMs were thawed and plated using the STEMdiff™ Cardiomyocyte Support Medium for 24 hours. After 24 hours, the spent medium was replaced with STEMdiff™ Cardiomyocyte Maintenance Medium. The medium was then replaced every 2 days.

FIGURE 5. hPSC-CMs produced with STEMdiff™ Cardiomyocyte Products are sensitive to HERG Potassium channel and L-type Calcium channel blockers



(A, B) Microelectrode Array (Maestro MEA, Axion Biosystems) recordings of hPSC-CMs (Day 27) show characteristic electrical profiles and drug response to E4031 (10 nM) and Nifedipine (300 nM). E4031 prolonged (A, gray line) and Nifedipine (B, gray line) shortened the repolarization, respectively.

Conclusions

- The STEMdiff™ Cardiomyocyte Products provide a robust and standardized complete workflow to produce and process hPSC-derived Cardiomyocytes.
- Easily and efficiently produce and maintain hPSC-derived Cardiomyocytes with the STEMdiff™ Cardiomyocyte Differentiation and Maintenance Kits. Confluent beating monolayers observed by day 15 and can be maintained up to a month or longer.
- Robust Cardiomyocyte differentiation from multiple PSC lines resulting in >80% cTnT-positive cells with a yield of >1 million cardiomyocytes from one well of a 12-well plate.
- Easy and fast dissociation of confluent beating monolayers of hPSC-derived cardiomyocytes in just 15 minutes using the STEMdiff™ Cardiomyocyte Dissociation Kit. Achieve >80% viable cardiomyocytes that can be used in downstream applications such as flow cytometry, cryopreservation, immunocytochemistry, or electrophysiology.
- Efficient cryopreservation and thawing of hPSC-derived Cardiomyocytes using the STEMdiff™ Freezing Cardiomyocyte Freezing Medium and STEMdiff™ Cardiomyocyte Support Medium.
- hPSC-derived Cardiomyocytes produced with the STEMdiff™ Cardiomyocyte Products show characteristic electrical waveforms and response to cardio-active drugs.