

# SSEA-4 PE EasySep® Positive Selection Kit for Selection of Pluripotent Stem Cells

Jennifer Antonchuk<sup>1</sup>, Alexandra Blak<sup>1</sup>, Jennifer Moody<sup>1</sup>, Andrea Eskander Afshari<sup>1</sup>, Mark Fairey<sup>1</sup>, Heather Drew<sup>1</sup>, Clive Glover<sup>1</sup>, Terry E. Thomas<sup>1</sup>, Allen C. Eaves<sup>1,2</sup> and Sharon A. Louis<sup>1</sup>

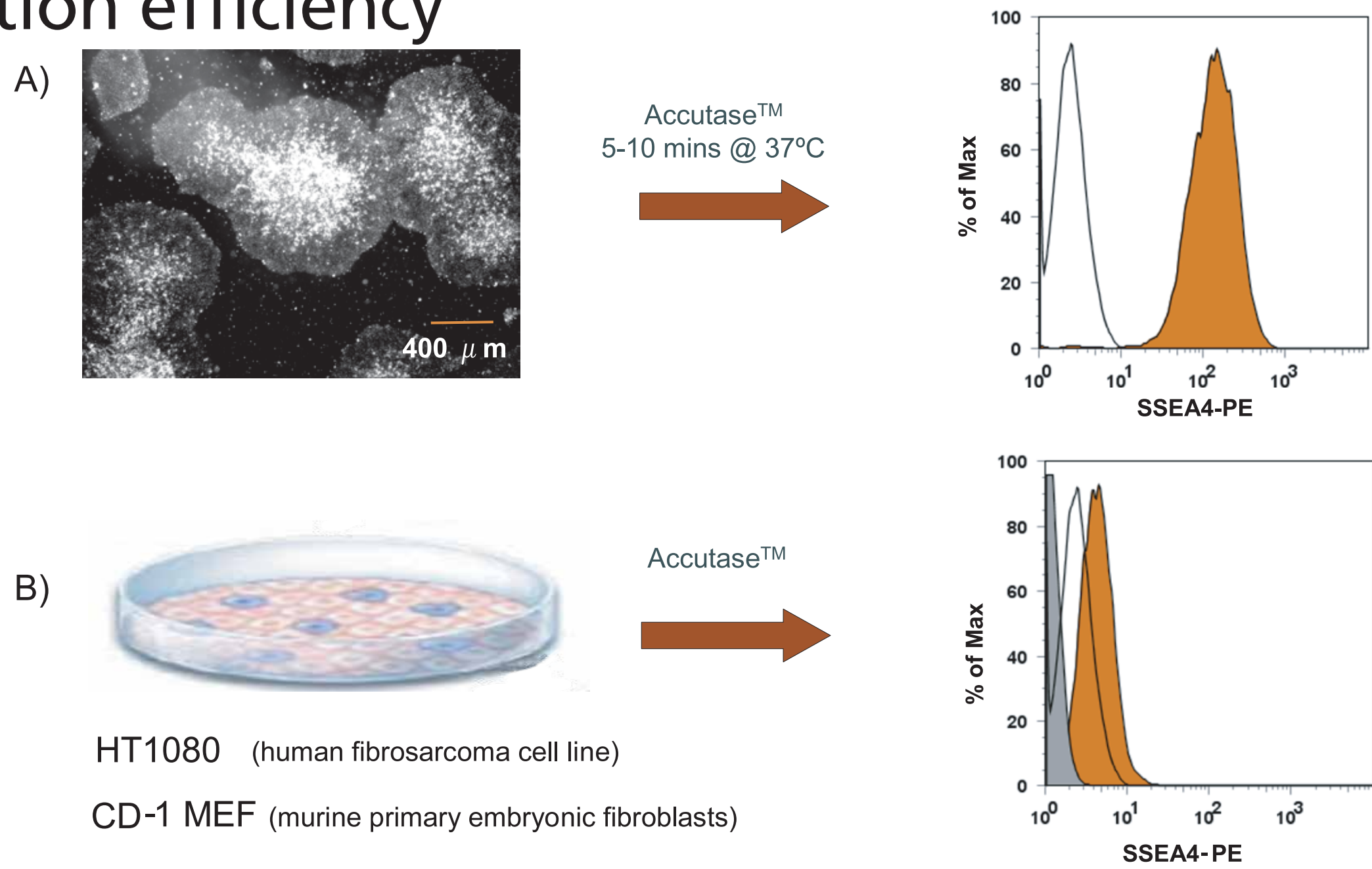
<sup>1</sup> STEMCELL Technologies Inc., Vancouver, BC, Canada  
<sup>2</sup> Terry Fox Laboratory, BC Cancer Agency, Vancouver, BC, Canada

## Abstract

Undifferentiated human pluripotent stem cells (PSCs) express SSEA-4 on their surface, and this expression is downregulated with differentiation, making it a reliable marker of pluripotency. We describe a rapid and simple method for the enrichment of PSCs by positive selection, that yields reliably high cell purities and recoveries. PSCs were isolated using immuno-magnetic, column-free positive selection (EasySep®), based on expression of SSEA-4. Importantly, the selected cells were able to reattach to Matrigel® coated plates in mTeSR®1 medium, and maintained characteristic undifferentiated morphology for at least 3 passages in mTeSR®1 following selection.

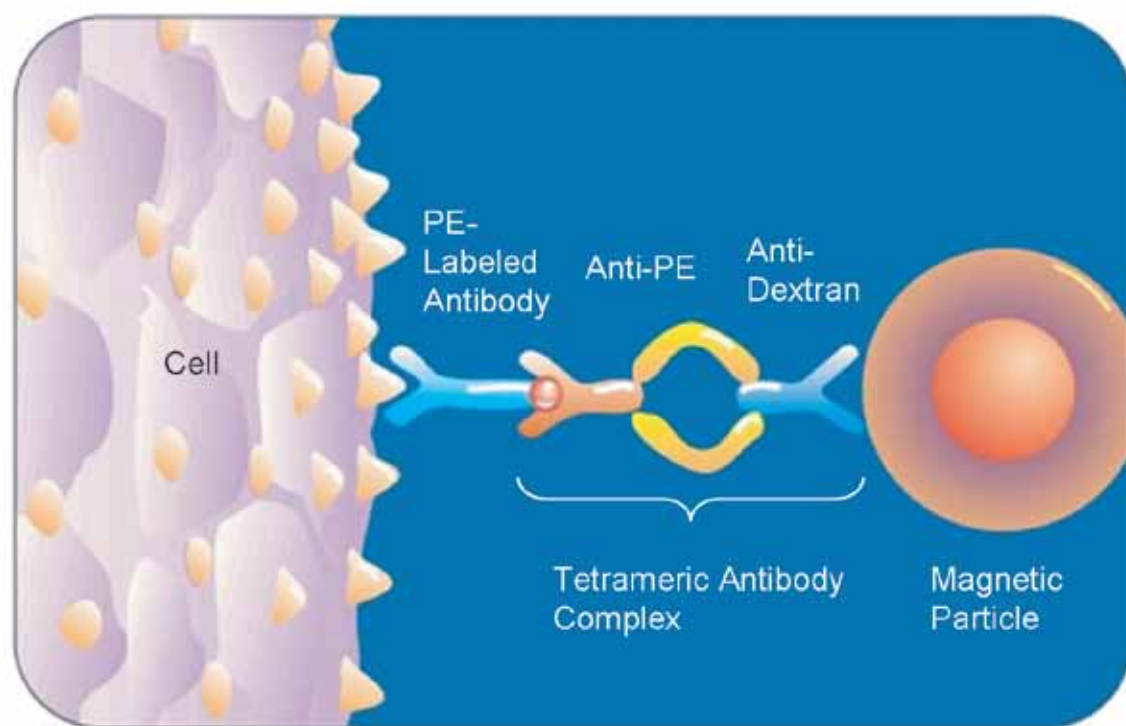
## Methods

Figure 1: SSEA-4<sup>+</sup> PSCs and SSEA-4<sup>-</sup> non-PSCs used in mixed populations to test selection efficiency



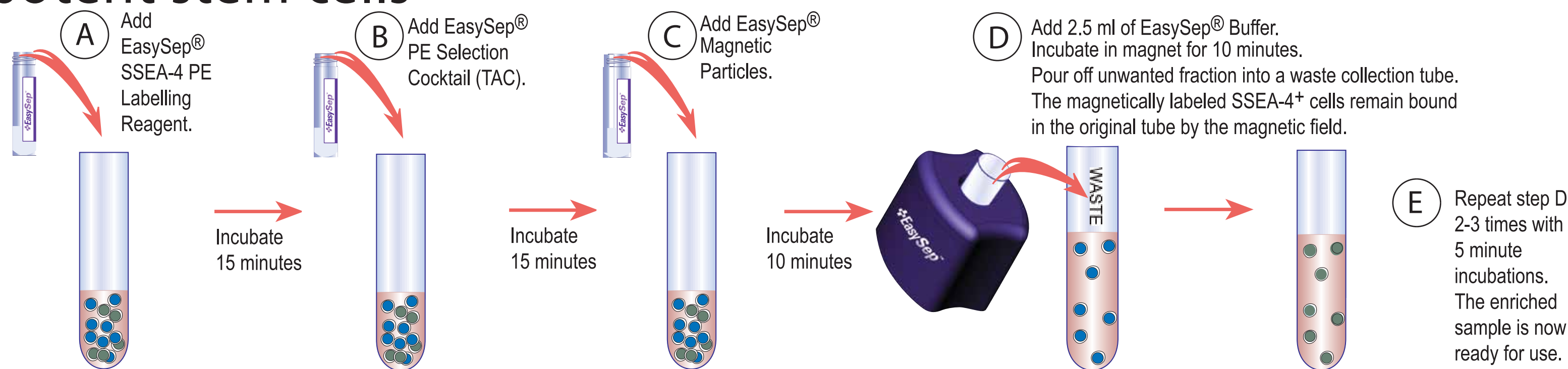
A) Single cell suspensions of H9 hES cells were obtained by incubating with Accutase®, and were confirmed by FACS analysis to be SSEA-4 positive (right: orange=H9 stained with SSEA-4 PE; white=isotype control). B) Single cell suspensions of non-PSC control cells: HT1080 fibrosarcoma cells or CD-1 murine embryonic fibroblasts (MEF) were obtained by incubating with Accutase® and were confirmed to be SSEA-4 negative (right: grey=MEF; orange=HT1080; white=isotype control).

Figure 2: EasySep® labeling of human pluripotent stem cells



SSEA-4 expressing cells are specifically labeled with anti-SSEA-4 PE antibody, followed by bispecific tetrameric antibody complex (TAC) and dextran-coated magnetic particles. Using the EasySep® magnet, the desired SSEA-4<sup>+</sup> cells can then be retained, and unwanted cells removed from the suspension.

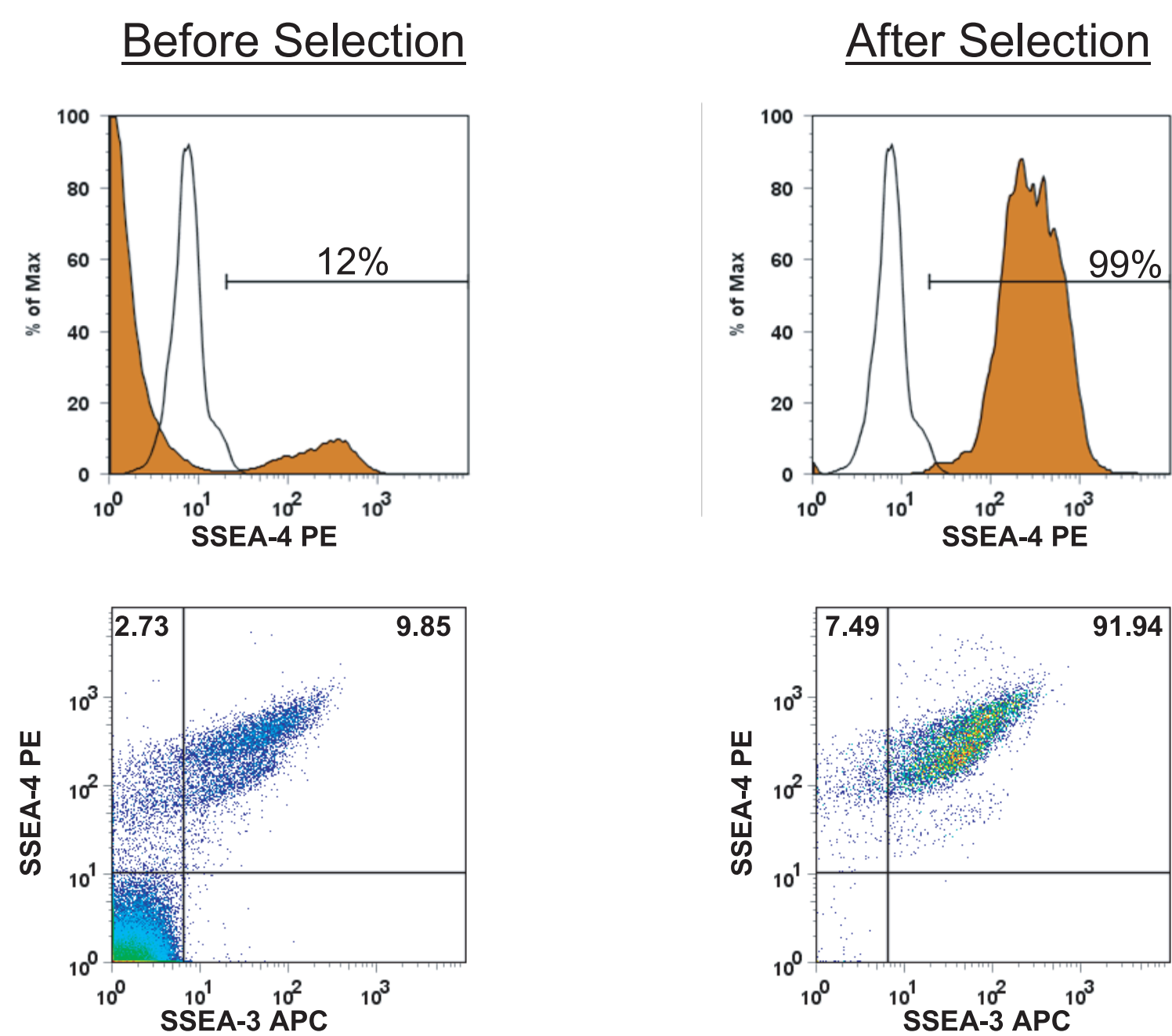
Figure 3: EasySep® procedure for column-free enrichment of human pluripotent stem cells



The EasySep® SSEA-4 PSC positive selection kit is designed to isolate human PSCs which express SSEA-4 (STEMCELL Cat #18165). (A-C) Single cell suspension from PSC-containing culture is incubated sequentially with EasySep® SSEA-4 PE Labelling Reagent, EasySep® PE Selection Cocktail (TAC), and EasySep® Magnetic Particles. D) After incubation in the magnet for 10 minutes, unlabelled cells are poured off into a waste collection tube while the magnetically labeled SSEA-4<sup>+</sup> cells are held in the original tube by the EasySep® magnet. E) 2-3 additional rounds of magnetic cell separation enable the isolation of highly purified PSCs that are ready to use in functional assays. The entire procedure requires only 60 minutes.

## Results

Figure 4: SSEA-4 EasySep® enrichment FACS profiles



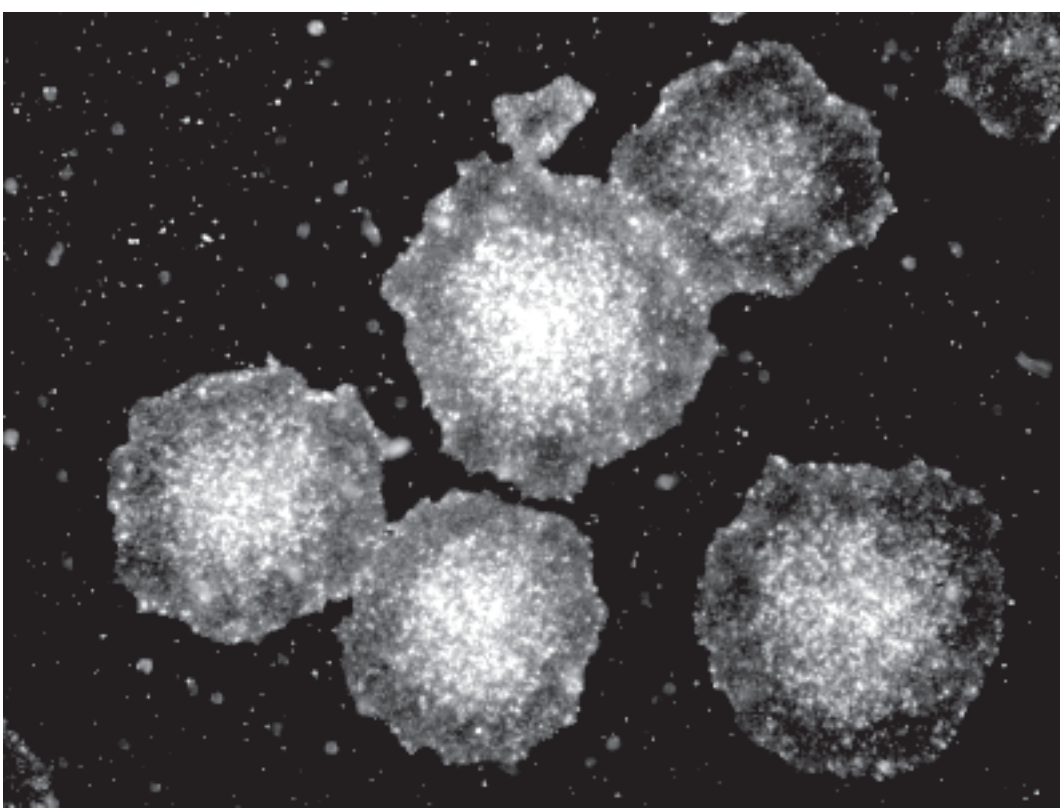
Cells were mixed at 1:9 ratio (10<sup>6</sup> PSCs and 9x10<sup>6</sup> non-PSCs), to simulate a mixed PSC population. Shown here is a representative experiment starting with a mixed population containing 12% SSEA-4<sup>+</sup> cells and resulting in 99% SSEA-4<sup>+</sup> cells after 4 rounds of separation (top: orange= SSEA-4 PE; white= isotype control, 1.24%<sup>+</sup>). Selected cells also co-stain with pluripotency marker SSEA-3 (bottom).

Table 1: Purity and recovery of SSEA-4<sup>+</sup> PSCs enriched by EasySep® positive selection using three or four rounds of selection

	# of Replicates	Purity (%)	Recovery (%)
3 rounds of selection	6	72 ± 6	63 ± 26
4 rounds of selection	7	88 ± 4	28 ± 15

Values expressed as mean ± SD  
Purity and recovery determined by flow cytometry. All samples gated on viable (PI negative) cells.

Figure 5: Selected PSCs maintain undifferentiated colony morphology for at least 3 passages after EasySep® selection



PSCs obtained after EasySep® selection from a mixed population of 10% hES cells and 90% MEF or HT1080 cells were plated onto Matrigel®-coated plates in mTeSR®1 medium. The cells were able to establish colonies with typical hES morphology, lacking overt signs of differentiation or noticeable fibroblast outgrowths.

## Conclusions

- PSCs can be rapidly isolated from mixed populations using column-free positive selection.
- High recovery and high purity can be achieved, by performing 3 or 4 rounds of EasySep® separation.
- Enriched PSCs can be used to re-seed PSC cultures to give normal colonies.



THE CELL EXPERTS™   WWW.STEMCELL.COM	
IN NORTH AMERICA	TOLL-FREE T. 1 800 667 0322 TOLL-FREE F. 1 800 567 2899 T. 1 604 877 0713 F. 1 604 877 0704 E. INFO@STEMCELL.COM TOLL-FREE T. 00 800 7836 2355 TOLL-FREE F. 00 800 7836 2300
IN EUROPE	T. +33 (0)4 76 04 75 30 F. +33 (0)4 76 18 99 63 E. INFO.EU@STEMCELL.COM
IN AUSTRALIA	TOLL-FREE T. 1 800 060 350 F. +61 (03) 9338 4320 E. INFO.AUS@STEMCELL.COM
IN SINGAPORE	T. (65) 6776 7754 F. (65) 6776 7114 E. INFO.SG@STEMCELL.COM