

A RAPID AND SIMPLE METHOD FOR THE ISOLATION OF HIGHLY FUNCTIONAL PLASMACYTOID DENDRITIC CELLS FROM HUMAN PERIPHERAL BLOOD

Andy I. Kokaji¹, Siobhan Holland¹, Victoria S. Stronge¹, Nicholas J. Megjugorac², Grant E. Gallagher², Maureen A. Fairhurst¹, Benoit Guilbault¹, Grant Gallagher², and Terry E. Thomas¹

¹ STEMCELL Technologies Inc., Vancouver, BC, Canada

² Genetic Immunology Laboratory, HUMIGEN LLC., Institute for Genetic Immunology, Hamilton, NJ, USA

ABSTRACT

Plasmacytoid DCs (pDCs) play an important role in antiviral immunity, secreting large amounts of type I and type III interferon (IFN) upon stimulation by live virus or toll-like receptor (TLR) agonists. pDCs comprise only a small fraction of total circulating peripheral blood mononuclear cells (PBMC) and current protocols for their isolation are time-consuming, involve multiple steps, and often require special equipment. We describe a rapid and simple method for the enrichment by negative selection of highly functional pDCs from normal blood that yield excellent cell purities and recoveries. Briefly, PBMCs are isolated by Ficoll-Paque PLUS density gradient sedimentation and pDCs are isolated using immuno-magnetic, column-free, negative selection (EasySep[®]). Our EasySep[®] technology involves specifically labeling unwanted cells with a cocktail of bispecific tetrameric antibody complexes (TAC) and dextran-coated magnetic particles. Using a magnet, the magnetically-labeled unwanted cells are removed from the unlabeled pDCs. Using our RoboSep[®] cell separator, the entire procedure can be fully automated. Flow cytometric assessment of EasySep[®]-isolated pDCs (Lin⁻, HLA-DR⁺, BDCA-4⁺) demonstrate purities of 93.8% ± 3.8% with cell recoveries of 65.6% ± 16.2% (n=9). Functional assessment of EasySep[®] and RoboSep[®]-isolated pDCs demonstrated robust secretion of type I and type III IFN following treatment with TLR agonists and live viruses.

METHODS

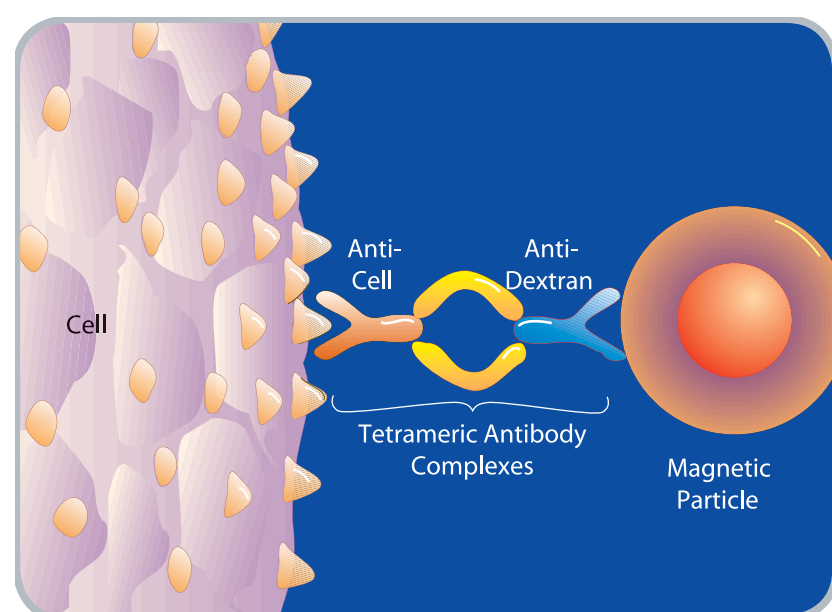
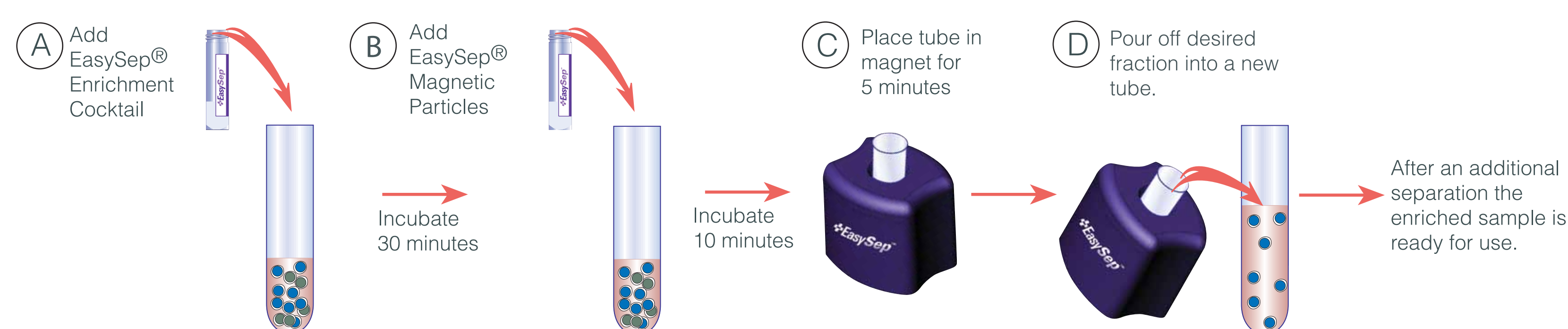


FIGURE 1: EasySep[®] labeling of human PBMC

Unwanted cells are specifically labeled with dextran-coated magnetic particles using bispecific TAC. These TAC recognize both dextran and the cell surface antigen expressed on the unwanted cells. Magnetically-labeled cells are then separated from unlabeled target cells using the EasySep[®] procedure.

FIGURE 2: EasySep[®] procedure for column-free enrichment of plasmacytoid dendritic cells



The EasySep[®] Human pDC Enrichment Kit is designed to isolate untouched pDCs from fresh or previously frozen PBMC samples by depletion of all other non-pDCs (STEMCELL Technologies, Inc. Cat #19062). A) PBMCs are incubated for 30 minutes with the EasySep[®] pDC enrichment cocktail; B) followed by a 10 minute incubation with EasySep[®] magnetic particles. C) The tube containing the labeled PBMCs is then placed in an EasySep[®] magnet for 5 minutes. D) EasySep[®] enriched pDCs are poured off into a new tube while the magnetically labeled non-pDCs cells are held in the original tube by the EasySep[®] magnet. An additional round of magnetic cell separation enables the isolation of highly purified pDCs that are ready to use in functional assays. The entire procedure requires less than 1 hour and can be fully automated with the RoboSep[®] cell separator.

RESULTS

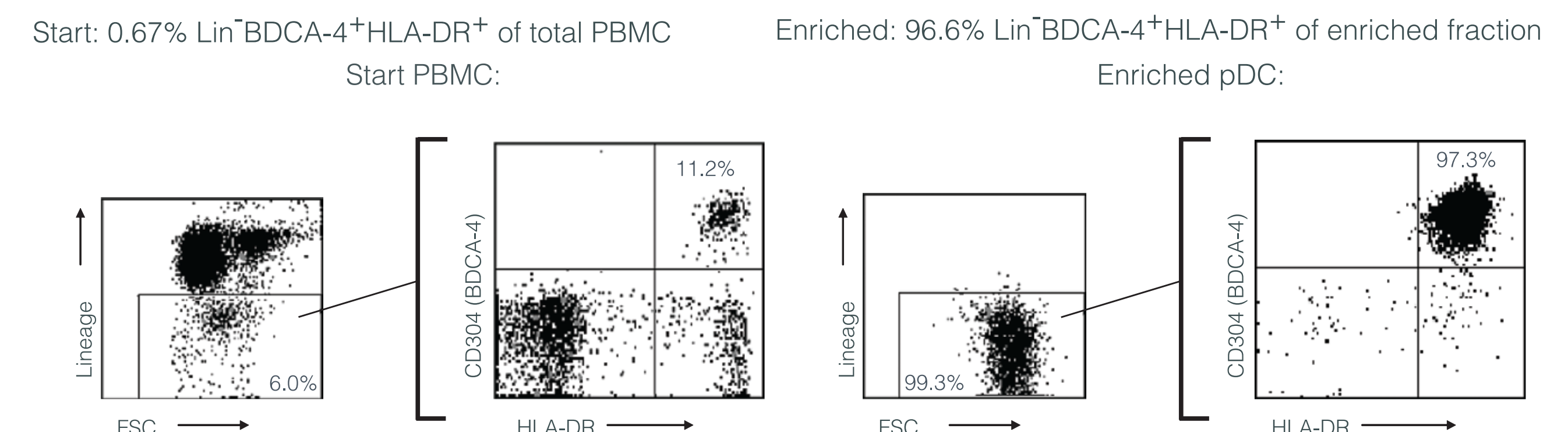
TABLE 1: Purity and recovery of Lin⁻HLA-DR⁺CD304⁺ pDCs enriched by negative selection from PBMC using EasySep[®] and RoboSep[®]

	n	Purity	Recovery
EasySep [®]	9	93.8 ± 3.8	65.6 ± 16.2
RoboSep [®]	5	92.8 ± 2.6	76.8 ± 20.6

Values expressed as means ± SD

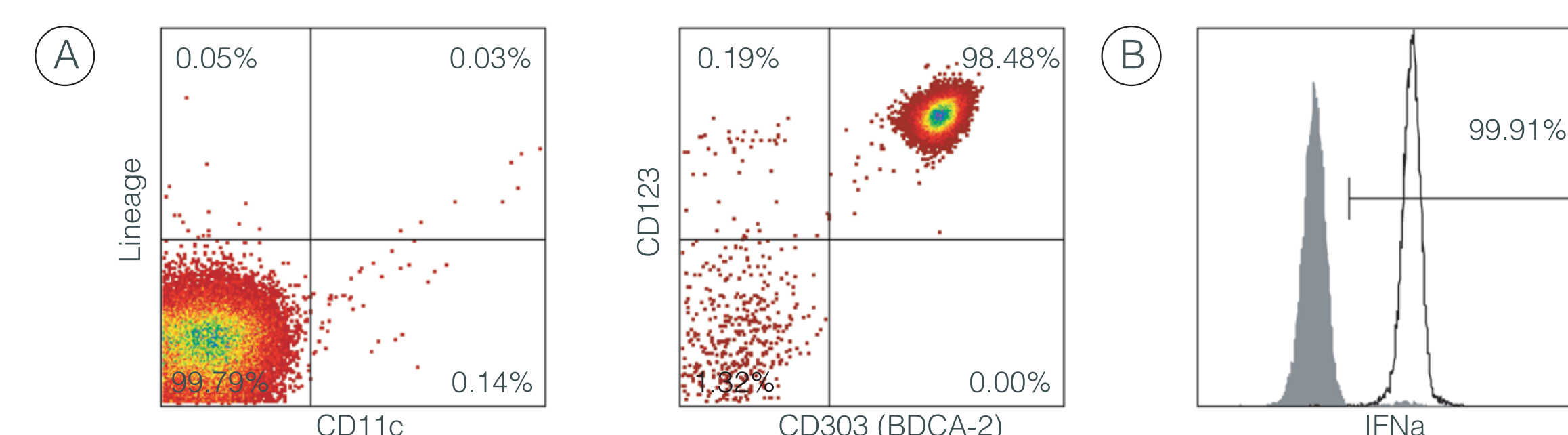
Purity and recovery determined by flow cytometry. All samples gated on viable (PI negative) cells.

FIGURE 3: Typical EasySep[®] and RoboSep[®] enrichment profiles



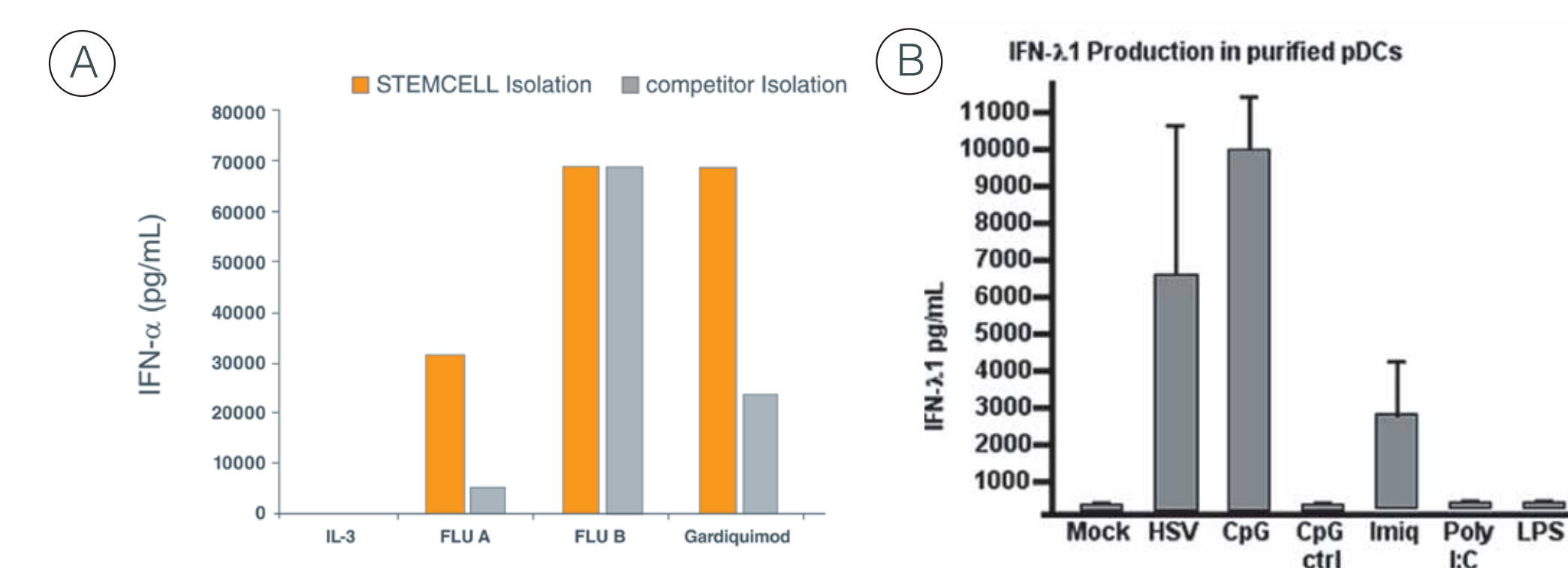
Starting with 0.2 - 0.9% pDCs in PBMC, the pDC content of the enriched fraction typically ranges from 87- 97% purity based on the pDC phenotype of lineage (CD3, CD14, CD16, CD19, CD20, CD34, CD56) negative, HLA-DR positive, and CD304 (BDCA-4) positive.

FIGURE 4: Virtually all EasySep[®] enriched pDCs express IFN α following *in vitro* TLR7 stimulation



Untouched human pDCs secrete IFN α following *in vitro* activation. A) EasySep[®]-enriched human pDCs were analyzed for the expression of lineage markers, CD11c, CD123 and CD303 (BDCA-2). B) EasySep[®]-enriched pDCs were cultured for 20 hours in serum-free media supplemented with 5 μ g/mL Imiquimod-R837, a small synthetic TLR7 ligand. Monensin was added during the final 2 hours of culture followed by flow cytometric assessment of intracellular IFN α . Open and shaded histograms represent staining with anti-human IFN α (clone: MMHA-11) or an isotype control, respectively.

FIGURE 5: EasySep[®] or RoboSep[®]- isolated pDCs secrete Type I and Type III IFN following stimulation with various stimuli



A) Compared to competitor isolated pDCs, RoboSep[®]-isolated pDCs secrete increased levels of IFN α following treatment with type A Influenza or a TLR7 agonist, Gardiquimod[™]. Data kindly provided by Dr. Michelle Gill from the Department of Pediatrics, University of Texas Southwestern Medical Center. B) EasySep[®]-isolated pDCs were incubated with various TLR agonists for 24 h and assayed for production of IFN λ 1 by ELISA.

CONCLUSIONS

- UNTOUCHED HUMAN PDCS CAN BE RAPIDLY ISOLATED FROM PBMC SAMPLES USING COLUMN-FREE NEGATIVE SELECTION.
- EXCELLENT PURITY AND RECOVERY CAN BE ACHIEVED WITH THE EASYSEP[®] AND ROBOSEP[®] PLATFORMS.
- VIRTUALLY ALL EASYSEP[®] ENRICHED PDCS EXPRESS IFN α FOLLOWING *IN VITRO* TLR7 STIMULATION WITH IMIQUIMOD-R837 IN SERUM-FREE MEDIA.