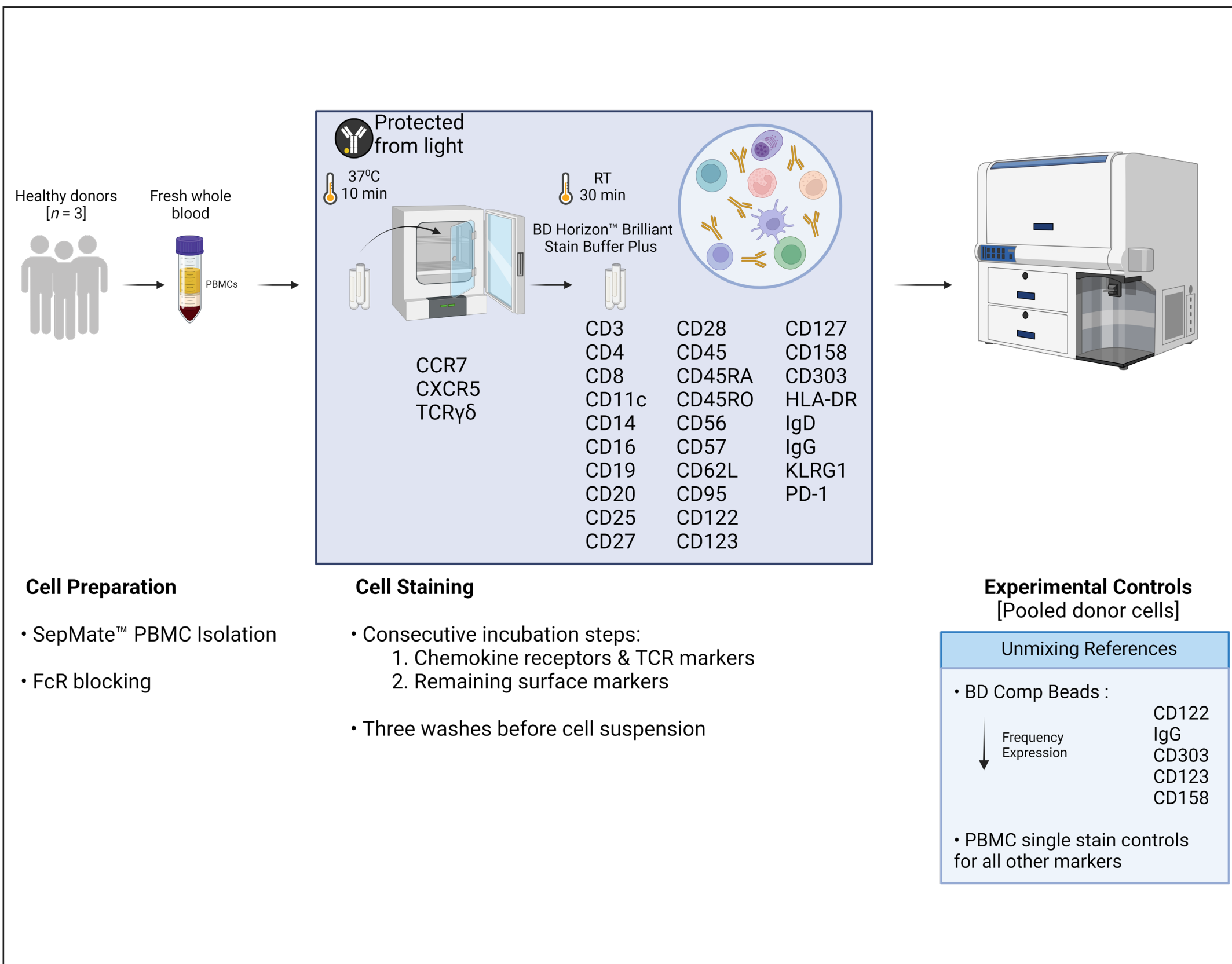


Abstract

In the rapidly advancing field of flow cytometry, cross standardization of flow cytometers remains a challenge. Cross standardization ensures consistency of data generated on separate cytometers. In this study, we demonstrate the transfer of application settings from the BD FACSymphony™ A5 SE Cell Analyzer to the BD FACSymphony™ S6 SE Cell Sorter, enabling convenient transfer of a 33-color panel between two spectral-enabled flow cytometers.

First, the BD FACSymphony™ A5 SE Cell Analyzer and the BD FACSymphony™ S6 SE Cell Sorter were checked for performance and characterized using the BD® Cytometer Setup and Tracking (CS&T) System. Second, both instruments were set up with optimized voltage values obtained from a volttration analysis. Third, Median Fluorescence Intensity (MFI) from mid-peak rainbow fluorescent particles was recorded for each of the 48 PMT detectors on the BD FACSymphony™ A5 SE Cell Analyzer. Next, new application settings were created on the BD FACSymphony™ S6 SE Cell Sorter with voltages set to match the MFI values from the BD FACSymphony™ A5 SE Cell Analyzer. Finally, CD4 samples were recorded on each instrument for confirmation of similar MFI values. Following cross standardization, the 33 -color immunophenotyping panel designed on the BD FACSymphony™ A5 SE Cell Analyzer was directly transferred to the matched BD FACSymphony™ S6 SE Cell Sorter, facilitating the physical isolation of target cell populations of interest.

Methods



Fresh human PBMCs were derived from whole blood obtained from healthy donors. PBMCs were isolated by Ficoll™ gradient centrifugation and resuspended in DPBS (1X). The staining solution comprised a cocktail of antibodies as per the specified test volumes (Table 1) suspended in BD Horizon™ Brilliant Stain Buffer Plus. Cells were stained by first pre-staining the freshly isolated PBMCs with antibodies against TCRγδ, CD185 and CD197 for 10 min at 37 °C. The antibody cocktail was then added to the pre-stained peripheral blood mononuclear cell (PBMCs) and the mixture was incubated for 30 min at room temperature in the dark. Single-stained controls were processed in parallel with the full panel. Cells were washed twice using cold BSA wash buffer (PBS and 0.01% Na₃). Cells were resuspended in wash buffer and stored on ice, protected from light, until acquisition. Viability dye 7-AAD was added at the specified test volume 5 min before acquisition on the flow cytometer. Optimized gain settings were derived for the BD FACSymphony™ A5 SE Cell Analyzer to generate Application Settings. Application Settings were then applied to maintain consistent fluorescence intensity values across experiments.

Results

1 Instrument configuration and reagent selection

Leukocyte	Laser	Marker	Fluorochrome	Catalog #	Clone	Compensation Control	Test size (μl or μg /test)
T cell	UV	CD45RA	BUV395	740315	5H9	PBMCs	0.06 μg
		Auto F	BUV496	749888	UCHL1	PBMCs	0.5 μg
		CD45RO	BUV563	748851	3G8	PBMCs	0.5 μg
		CD16	BUV661	741559	RF8B2	PBMCs	0.06 μg
		CD185/CXCR5	BUV737	612815	CD28.2	PBMCs	0.25 μg
		CD127	BUV805	748486	HIL-7R-M21	PBMCs	0.5 μg
		CD25	BV421	567485	BC96	PBMCs	2.5 μl
		CD20	V450	561164	L27	PBMCs	2.5 μl
		CD62L	BV480	566174	DREG-56	PBMCs	0.125 μg
		CD122/IL2-RB	BV510	747741	Mik-β3	Beads	2.5 μl
B cell	Violet	IgG	BV605	563246	G18-145	Beads	2.5 μl
		CD56	BV650	565057	NCAM16.2	PBMCs	0.125 μg
		CD303	BV711	748002	V24-785	Beads	1.25 μg
		KLRG1	BV750	753692	Z7-205:mAb	PBMCs	1.25 μg
		CD123	BV786	564196	7G3	Beads	0.125 μg
		CD57	FITC	555619	NK-1	PBMCs	0.03 μg
		CD4	RB545	569183	SK3	PBMCs	0.06 μg
		CD11c	BB630-P2	Custom	B-LY6	PBMCs	0.5 μg
		CD27	BB660-P2	Custom	M-T271	PBMCs	0.0125 μg
		CD45	PerCP	340665	2D1	PBMCs	10 μl
NK cell	Blue	CD279/PD-1	BB700	566460	EH12.1	PBMCs	2.5 μl
		CD3	BB755-P	Custom	UCHT1	PBMCs	0.025 μg
		CCR7/CD197	RB780	568748	2-LT-A	PBMCs	1 μg
		CD19	PE	555413	HIB19	PBMCs	5 μl
		CD158	RV586	753232	HP-MA4	Beads	0.125 μg
		CD14	PE-CF594	562335	MqP99	PBMCs	0.125 μg
		CD95	PE-Cy5	559773	DX2	PBMCs	20 μl
		7-AAD	TCRγδ	559925	11F2	PBMCs (10% heat killed)	5 μl
		HLA-DR	APC	559868	TU36	PBMCs	1.25 μg
		IgD	R718	567993	IA6-2	PBMCs	20 μl
Dendritic cell	Yellow Green	CD3	BB700	566460	UCHT1	PBMCs	0.025 μg
		CD19	PE	555413	HIB19	PBMCs	5 μl
		CD158	RV586	753232	HP-MA4	Beads	0.125 μg
		CD14	PE-CF594	562335	MqP99	PBMCs	0.125 μg
		CD95	PE-Cy5	559773	DX2	PBMCs	20 μl
		7-AAD	TCRγδ	559925	11F2	PBMCs (10% heat killed)	5 μl
		HLA-DR	APC	559868	TU36	PBMCs	1.25 μg
		IgD	R718	567993	IA6-2	PBMCs	20 μl
		CD8	APC-H7	641409	SK1	PBMCs	0.03 μg
		Monocyte	Red	CD3	BB700	566460	UCHT1
CD19	PE			555413	HIB19	PBMCs	5 μl
CD158	RV586			753232	HP-MA4	Beads	0.125 μg
CD14	PE-CF594			562335	MqP99	PBMCs	0.125 μg
CD95	PE-Cy5			559773	DX2	PBMCs	20 μl
7-AAD	TCRγδ			559925	11F2	PBMCs (10% heat killed)	5 μl
HLA-DR	APC			559868	TU36	PBMCs	1.25 μg
IgD	R718			567993	IA6-2	PBMCs	20 μl
CD8	APC-H7			641409	SK1	PBMCs	0.03 μg
CD14	PE-CF594			562335	MqP99	PBMCs	0.125 μg

Table 1. Instrument configuration and reagent selection. This 33-color panel allows identification and sorting of specific human (PBMC) populations. Careful evaluation of single-color controls identified superior performance in spectral unmixing using BD® CompBeads Particles in place of cell based controls for five colors (BV510, BV605, BV711, BV786 and RY586) in this panel.

2 Identification of classical human immune cell subsets circulating in human blood

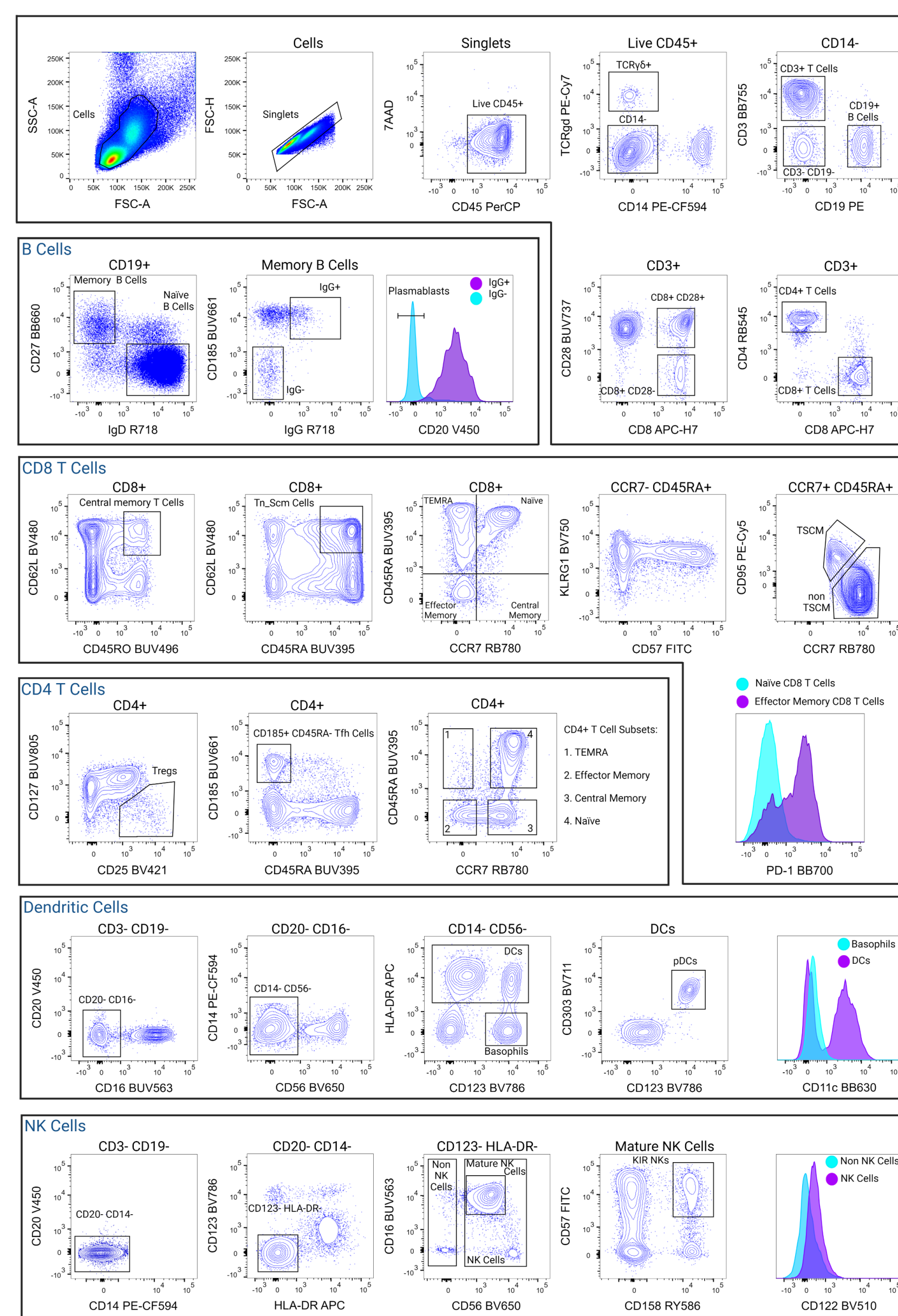
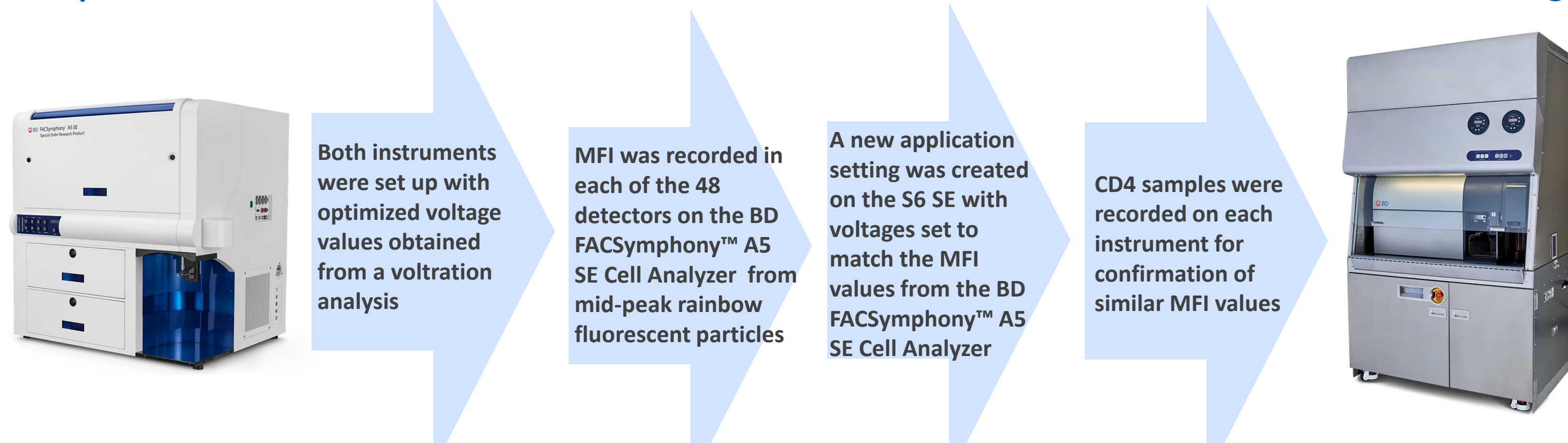


Figure 2. T cells, B cells, DCs and NK cells were gated on CD45+ cells by excluding 7AAD+ inviable cells and doublets. B cells were identified as CD19+ followed by identification of Naïve/Mature CD27-IgD+ B cells. Plasmablasts (CD27+ CD20-) were identified by assessing IgD- IgG- CD185- B cells. Classical T cells were identified as either CD4+, CD8+ or TCRγδ+, followed by identification of well-characterized T cell subsets based on the expression of CD62L and CD45RA or CD45RO (central memory and TN/SCM naïve and stem memory T cells), CD45RA and CCR7 (naïve, central memory, effector memory and terminally differentiated effector memory cells), CD95 and CCR7 (stem memory T cells TSCM), CD127 and CD25 (Tregs), and CD185 and CD45RA (T follicular helper cells). PD1 expression was assessed in CD8+ naïve and effector memory cells, and KLRG1 expression was assessed in CD8+ TEMRA cells. Classical DCs were identified as CD3- CD19- CD20- CD16- CD14- CD56- HLA-DR+, followed by identification of the pDC subset exclusively as CD303+CD123+. Basophils were identified as CD3- CD19- CD20- CD16- CD14- CD56- HLA-DR- CD123+. NK cells were identified as CD3- CD19- CD20- CD14- CD123- HLA-DR-. Mature and immature NK cells were distinguished based on the expression of CD16 and CD56 followed by identification of KIR-NK cells as CD57+CD158+ mature NK cells. NK cells and non-NK cells were assessed for the expression of CD122.

Results

3 High-dimensional UMAP analysis of multiparameter flow cytometry data for identification of donor variable cell subsets for cell sorting

4 Instrument Standardization



5 Transfer of the 33 color panel from the BD FACSymphony™ A5 SE Cell Analyzer to the BD FACSymphony™ S6 SE Cell Sorter for cell isolation

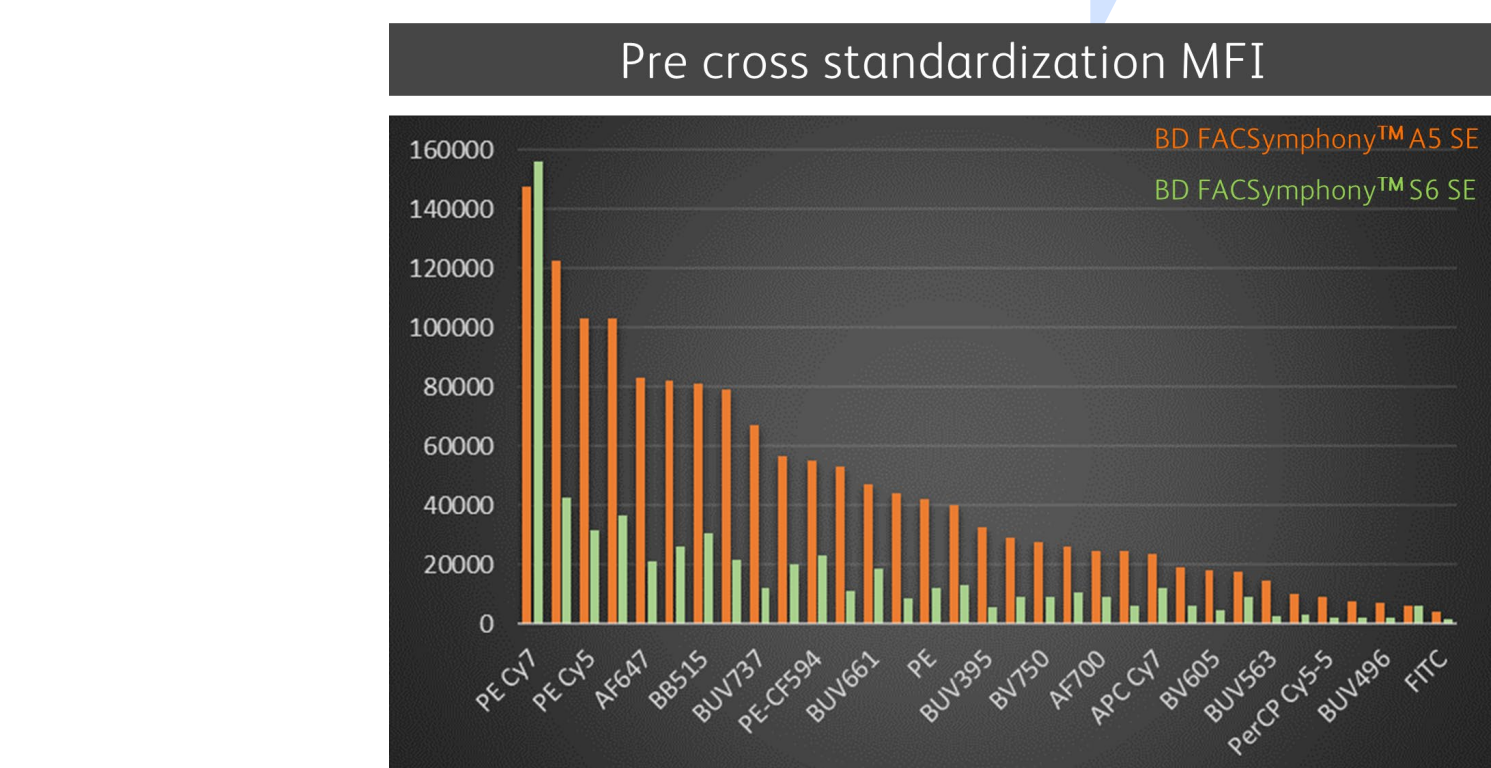
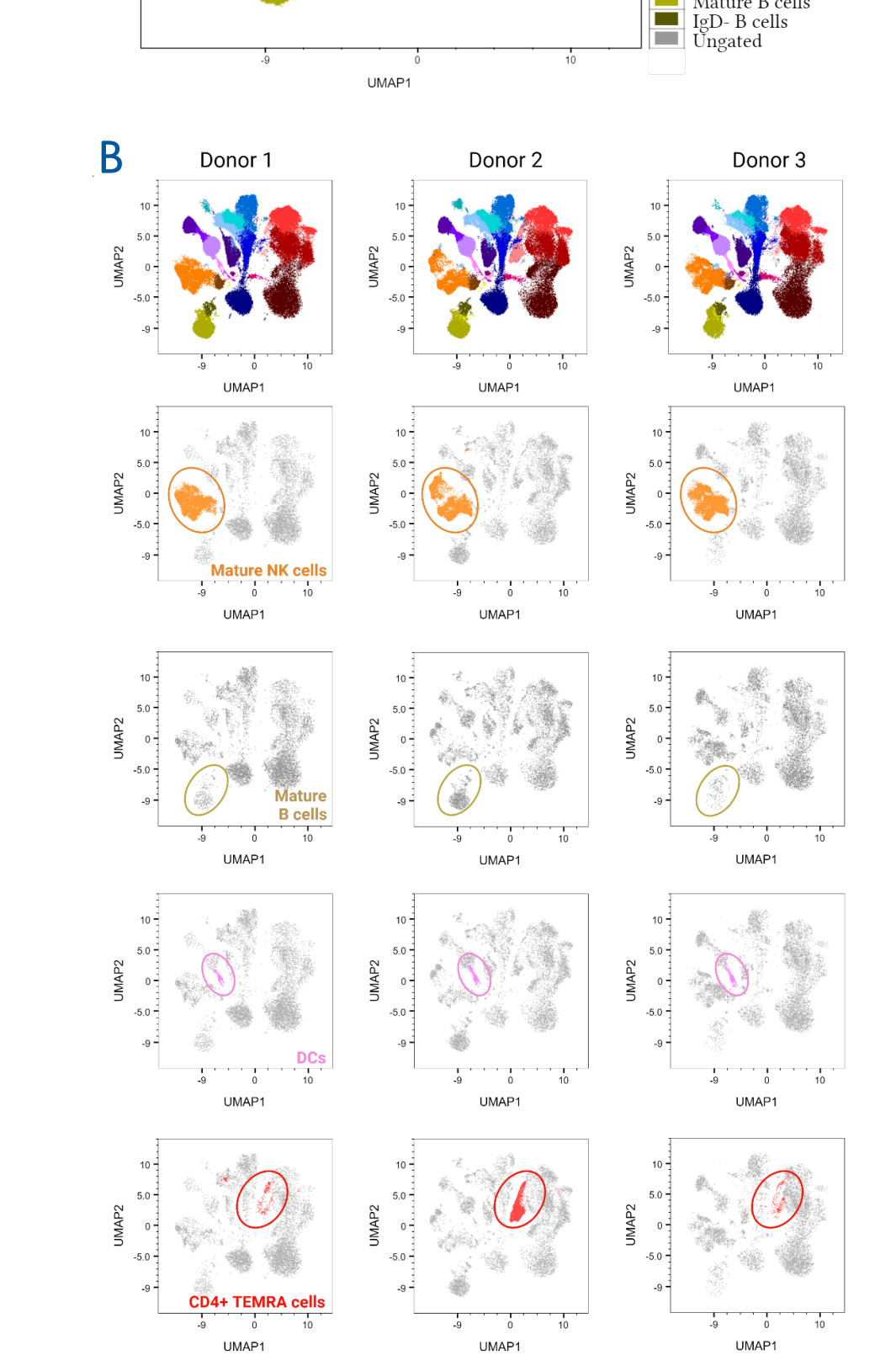
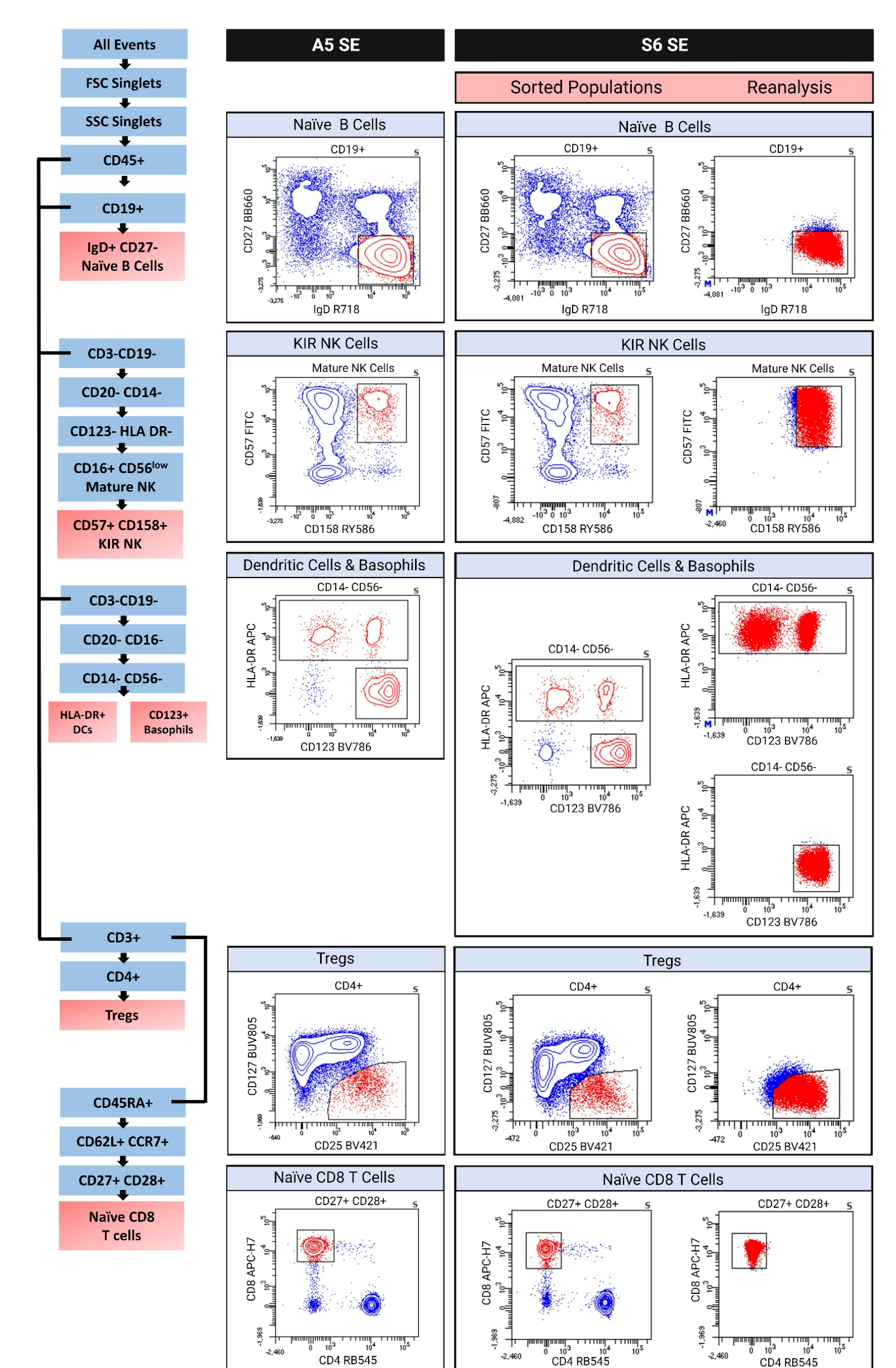


Figure 3A. Algorithmic analysis of multiparameter flow cytometry data was conducted using dimensionality reduction with UMAP and unsupervised clustering with FlowSOM* on concatenated CD45+ cells from n=3 donors. Clusters were annotated using the ClusterExplorer and Marker Enrichment Model tools and thereafter color-coded.

Figure 3B: UMAPs highlighted variations in the frequency of specific cell populations among the three donors.

Figure 4. MFI from CD4 samples recorded on the BD FACSymphony™ A5 SE Cell Analyzer and the BD FACSymphony™ S6 SE Cell Sorter prior to (left) and post (right) cross standardization.

Figure 5. Following cross-standardization, the panel was transferred to the BD FACSymphony™ S6 SE Cell Sorter. Representative data from the same sample demonstrate the comparable phenotypes of cell populations analyzed on the BD FACSymphony™ S6 SE Cell Sorter. Reanalysis of the sorted target populations, on the BD FACSymphony™ S6 SE Cell Sorter, reveals fidelity to the sort regions as gated on the parent population.

Conclusions

- Panel Design: We described a 33-color spectral panel that provides resolution of specific T, B, NK and DC cell populations in human healthy PBMC donor samples.
- Data analysis: High-dimensional UMAP analysis of multiparameter flow cytometry data allowed identification of donor variable cell subsets.
- Spectral unmixing: Live spectral unmixing of multiparameter flow cytometry panel allowed identification of cell subsets for cell isolation.

Class 1 Laser Product. For Research Use Only. Not for use in diagnostic or therapeutic procedures.

BD, the BD logo and [enter all product brand names in poster] are the trademarks of Becton, Dickinson or its affiliates. © 2024 BD. All rights reserved. BD-xxxxxx (v1.0) 0524