



High Purity Cell Isolation for Chimerism Analysis

Fully Automated, High Purity Cell Isolation for Chimerism Labs

Chimerism analysis is crucial to monitoring post-transplant outcome and determining appropriate post-transplantation therapy. Chimerism tests require **sensitive and reliable techniques** to detect the presence of donor leukocytes in the hosts' peripheral blood or bone marrow. Ensuring that cross-contamination between samples does not occur is critical to the success of downstream chimerism analyses.

RoboSep™, the **fully-automated** cell separator from STEMCELL Technologies Inc., **eliminates the risk of cross-contamination** during cell separations by using individual filter tips to process each cell sample. Based on **EasySep™**, the column-free immunomagnetic cell separation platform, RoboSep™ isolates **highly purified** cells from virtually any sample **in as little as 45 minutes**. Large routine laboratories worldwide rely upon RoboSep™ for high purity cell separation that enables clear and reliable downstream chimerism analysis.



www.stemcell.com/robosep

Ready·Sep·Go with RoboSep™

HIGH PURITY.

Achieve purities of up to 99% in as little as 45 minutes with minimal hands-on time.

VERSATILE.

Customize protocols for any cell type and sample source.

EFFICIENT.

Isolate multiple cell subsets from a single, low-volume sample.

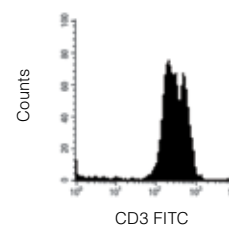
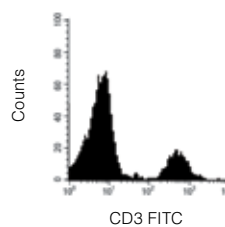
NO CROSS-CONTAMINATION.

Maintain sample integrity: Samples are processed with single-use disposable tips.

FACS Histogram Results with RoboSep™ HLA Buffy Coat / Whole Blood CD3⁺ Cell Selection Kit

Start*: 16% CD3⁺ Cells

Selected: 99% CD3⁺ Cells



*Red blood cells were removed by lysis prior to flow cytometry.



Scientists Helping Scientists™ | WWW.STEMCELL.COM

CATALOG #29173 | VERSION 2.0.1 | OCT 2011

TOLL-FREE T. 1 800 667 0322 • T. +1 604 877 0713 • TOLL-FREE F. 1 800 567 2899 • F. +1 604 877 0704
ORDERS@STEMCELL.COM • INFO@STEMCELL.COM • FOR FULL CONTACT DETAILS WORLDWIDE VISIT OUR WEBSITE

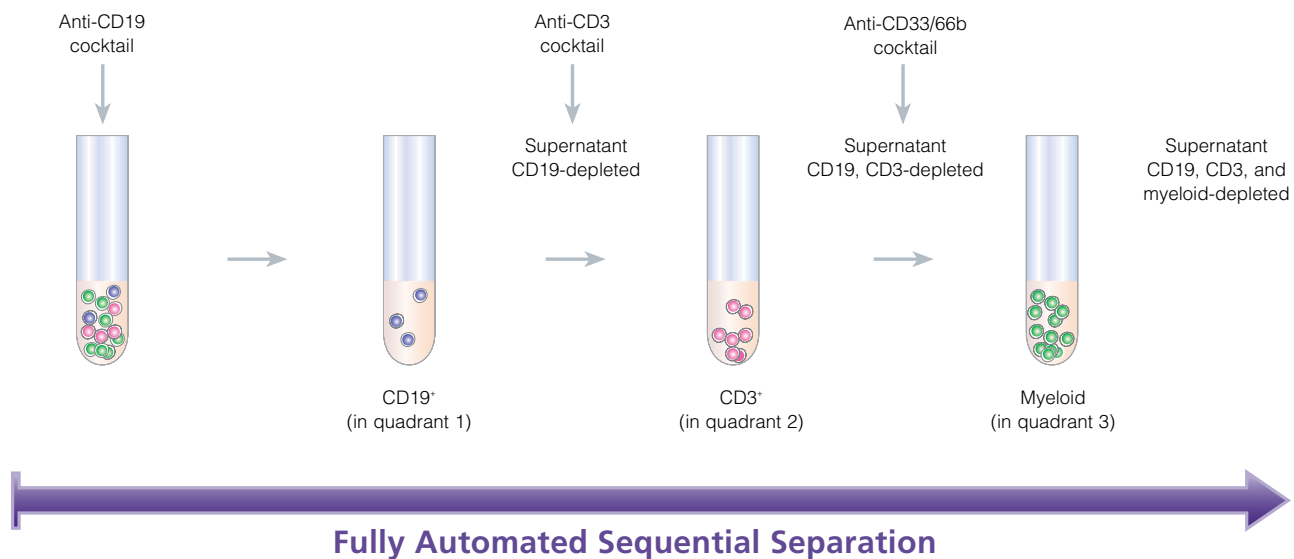
FOR RESEARCH USE ONLY. NOT FOR THERAPEUTIC OR DIAGNOSTIC USE.

Application

Fully Automated Separation of Multiple Cell Types from a Single Sample

Because chimerism analysis is typically performed on small blood samples, analysis of purified cell subsets often requires isolation of more than one cell type from a single starting sample. Sequential separation protocols from RoboSep™ can isolate up to four highly-enriched cell types from a single, undivided blood sample. Cells are recovered with high purity and in sufficient quantities for downstream flow cytometry and DNA analyses.

FIGURE 1. RoboSep™ sequential separation of B cells, T cells, and myeloid cells from a single sample.



Product Listing

CELL TYPE	CELL TYPE MARKER	RoboSep™	
		BUFFY COAT/WHOLE BLOOD ¹	PBMC, BM, SPLEEN, LN ²
Total Lymphocytes	CD3/CD19	18684HLARF	-
T Cells	CD3	18081HLARF	18051HLARF
B Cells	CD19 CD19/CD20	18084RF 18184HLARF	18054RF 18454HLARF
Myeloid Cells	CD15 CD33 CD33/66b	18681HLARF 18287HLARF 18683HLARF	- 18257RF -
Granulocytes	CD66b	18682RF	-
Monocytes	CD14	18088RF	18058RF
NK Cells	CD56	18085HLARF	18055RF
Hematopoietic Progenitor Cells	CD34	18086RF	18056RF

1. Kit also works on other red blood cell containing samples (i.e. cord blood, buffy coat, bone marrow).
2. PBMC = Peripheral Blood Mononuclear Cell; BM = Bone Marrow; LN = Lymph Node