

Multi-Color Flow Cytometry Panels for Immunophenotyping of Expanded T Cells

Introduction

Bio-Techne offers many immune cell therapy solutions, including fluorescent-conjugated antibodies for flow cytometry to assess T cell memory phenotype, T cell activation, and T cell exhaustion. Together with our Fluorokines™ fluorescent-labeled proteins, these panels will streamline the path to immune cell analysis during CAR-T cell development.

Immune cell therapy encompasses several approaches to harness the host immune response for treating immune disorders and cancer, including chimeric antigen receptor (CAR) T cell therapy, regulatory T cell (Treg) therapy and tumor infiltrating lymphocytes (TILs) therapy. Production of any therapeutic involving biological materials has a high level of inherent variability. Personalized cell therapies like CAR-T cell therapy add additional layers of complexity.

CAR-T cell therapy manufacturing is a multi-stage process that involves isolating T cells from donors and engineering them to express a chimeric antigen receptor (CAR) to target specific antigens on a patient's tumor cells. After expansion, these CAR-T cells are transfused back into the patient, resulting in T cell activation and destruction of the tumor cells. Throughout the process, it is critical to monitor the memory phenotype, activation status, and exhaustion markers on the T cells. The most common way to monitor T cell phenotype is by flow cytometry to ensure successful activation, expansion, CAR-T expression and functionality of the CAR-T cells post-infusion.

Our experts have selected antibody-fluorochrome combinations that work together for these flow cytometry panels to save you time and money developing a flow cytometry panel from scratch. In this application note, we'll show how Bio-Techne fluorescent-conjugated antibodies and Fluorokines™ can be used to monitor the frequency and phenotype of peripheral blood mononuclear cells (PBMCs) and expanded T cells in immune cell therapy workflows.

When to Use Multi-Color T Cell Panels in Immune Cell Therapy Workflow

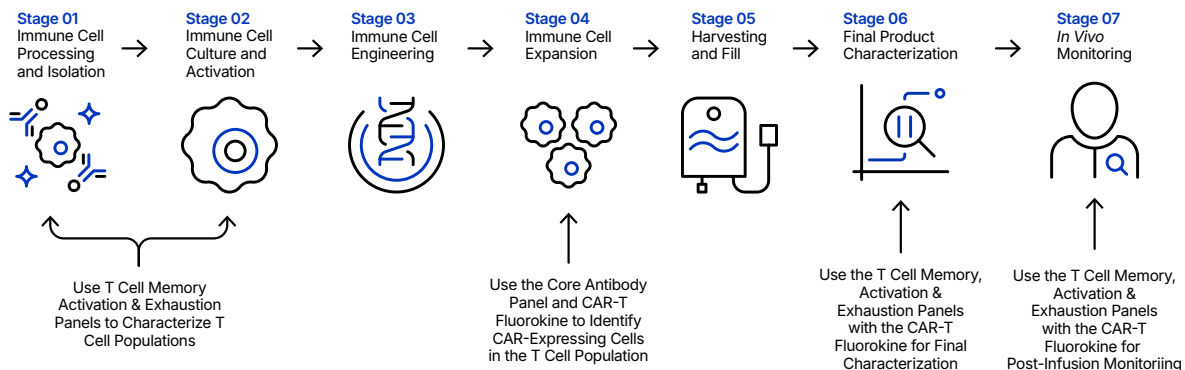


Figure 1: Overview of Patient-Specific Immunotherapy Workflow. Donor PBMC sample is processed for T cell selection (Stage 1) and enrichment (Stage 2). T cells are modified for chimeric antigen receptor (CAR) expression (Stage 3) to recognize tumor cell receptors and expanded in cell culture (Stage 4) before harvesting (Stage 5). The final CAR-T cell product is characterized (Stage 6) prior to re-introduction back into the patient (Stage 7).

Benefits of Using Multi-Color T Cell Panels

- Eliminate the guesswork from flow cytometry panel design by entrusting it to our team of antibody specialists.
- Enhance reproducibility and standardize immunophenotyping in your lab.
- Achieve reliable results with Bio-Techne's flow-validated antibody panels, featuring our rigorously tested R&D Systems™ antibodies.
- Compatible with Alexa Fluor® 488-conjugated Fluorokines and Near IR viability dyes.

Flow Cytometry Panel for Immunophenotyping T Cell Memory Markers

The frequency of various memory T cell subsets in a sample can significantly influence the expansion potential and functionality of the host immune response. Utilize this validated multi-color flow cytometry panel to phenotype human PBMCs for memory T cell markers, such as CD45RO and CCR7.

Flow Cytometry Panel for Immunophenotyping T Cell Activation Markers

T cells are activated when exposed to their specific antigen or external stimuli *in vitro*. Analyzing the expression patterns of early and late activation markers can enhance the understanding of the host immune response. Use this validated multi-color flow cytometry panel to identify activation markers, such as CD25, CD69, and CD38 on your T cells.

Flow Cytometry Panel for Immunophenotyping T Cell Exhaustion Markers

The expression of inhibitory receptors is essential for regulating immune responses. However, abnormal and prolonged expression of these receptors can cause immune dysfunction. Utilize this validated multi-color flow cytometry panel to analyze your T cells for exhaustion markers, such as PD-1, TIM-3, and LAG-3.

TABLE // 04

Selection of Fluorokines for CAR-Detection

| | | | | |
|----------------------------|----------------------|---------------|----------------|---------------|
| BCMA | CD19 | CD30/TNFRSF8 | CD40/TNFRSF5 | CD300e/LMIR6 |
| DLL3 | EGFR | EMMPRIN/CD147 | EpCAM/TROP1 | ErbB2/Her2 |
| Fc gamma RIIA/CD32a (R167) | Fc gamma RIIIA/CD16a | Glypican 3 | IL-13 R alpha2 | Mesothelin |
| MUC-1 | OX40 | PD-1 | Siglec-2/CD22 | Siglec-3/CD33 |
| TRAIL R2 | TSLPR | VEGFR2/KDR | | |

TABLE // 01

| Marker | Fluorochrome | Catalog # |
|--------|--------------------|----------------|
| CD3 | mFluor™ Violet 450 | FAB100MFV450 |
| CD4 | mFluor™ Violet 500 | FAB3791MFV500 |
| CD8 | Alexa Fluor® 700 | FAB1509N |
| CD45RA | mFluor™ Violet 610 | FAB11444MFV610 |
| CD45RO | Alexa Fluor® 647 | FAB10642R |
| CCR7 | PE | FAB197P |

TABLE // 02

| Marker | Fluorochrome | Catalog # |
|--------|--------------------|---------------|
| CD3 | mFluor™ Violet 450 | FAB100MFV450 |
| CD4 | mFluor™ Violet 500 | FAB3791MFV500 |
| CD8 | Alexa Fluor® 700 | FAB1509N |
| CD25 | mFluor™ Violet 610 | FAB1020MFV610 |
| CD38 | Alexa Fluor® 647 | FAB2404R |
| CD69 | PE | FAB23591P |

TABLE // 03

| Marker | Fluorochrome | Catalog # |
|--------|--------------------|----------------|
| CD3 | mFluor™ Violet 450 | FAB100MFV450 |
| CD4 | mFluor™ Violet 500 | FAB3791MFV500 |
| CD8 | Alexa Fluor® 700 | FAB1509N |
| LAG-3 | mFluor™ Violet 610 | FAB23193MFV610 |
| PD-1 | Alexa Fluor® 647 | FAB10863R |
| TIM-3 | PE | FAB2365P |

mFluor™ is used under license from AAT Bioquest, Inc. Alexa Fluor® is a registered trademark of Molecular Probes, Inc.

Flow Cytometry Panels Can Be Used to Detect T Cell Memory, Exhaustion, and Activation Markers in Human PBMCs

Initially, PBMCs are isolated from patient donors. Since the final output of CAR-T cells can be affected by the starting raw materials, the frequency of memory T cells, activated T cells and exhausted T cells by flow cytometry sets a baseline for the heterogeneity of the immune cell populations and, if isolating T cells before engineering, the overall purity of the sample (Figure 1: Stage 1).

Figure 2 shows the subpopulations of T cells expressing markers of memory (CD45RO/CD45RA), activation (CD69/CD38), and exhaustion (TIM-3/LAG-3) from a donor's PBMCs at Day 0, or Stage 1, of the immune cell therapy process.

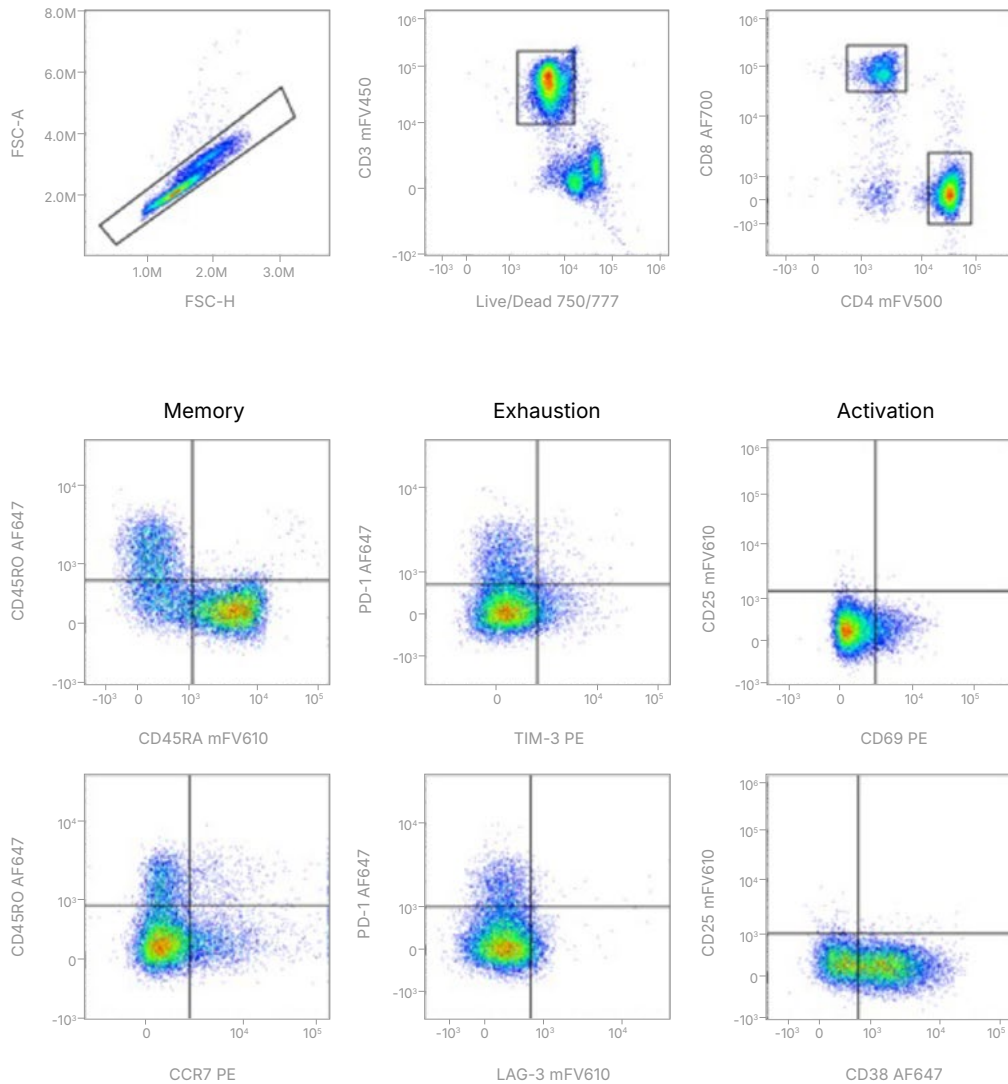


Figure 2: Characterization of PBMC Cells for T Cell Memory, Activation and Exhaustion with Multi-Color T Cell Panels using Flow Cytometry. Cryopreserved human (Day 0) PBMCs were thawed and stained with three flow cytometry panels to assess their naïve, memory, exhaustion, and activation phenotypes prior to T cell activation. PBMCs were stained with anti-human CD3 mFluor™ Violet 450, CD4 mFluor™ Violet 500, CD8 Alexa Fluor® 700, and Live/Dead 750/777 viability dye, followed by staining with three different panels to assess T cell 1) Memory (CD45RO Alexa Fluor® 647, CD45RA mFluor™ Violet 610, and CCR7 PE), 2) Exhaustion (PD-1 Alexa Fluor® 647, TIM-3 PE, and LAG-3 mFluor™ Violet 610), and 3) Activation (CD25 mFluor™ Violet 610, CD69 PE, and CD38 Alexa Fluor® 647).

Flow Cytometry Panels Identify Changes in Human T Cell Phenotypes After Activation

During T cell activation and cell culture expansion (Figure 1: Stage 2), immunophenotyping with flow panels is used to evaluate changes in T cell activation and overall changes in the distribution of T cell subtypes. These can be critical for understanding

the success of T cell engineering and output of the final product. Figure 3 shows representative staining using Memory T Cell, T Cell Exhaustion, and T Cell Activation flow cytometry panels from Bio-Techne.

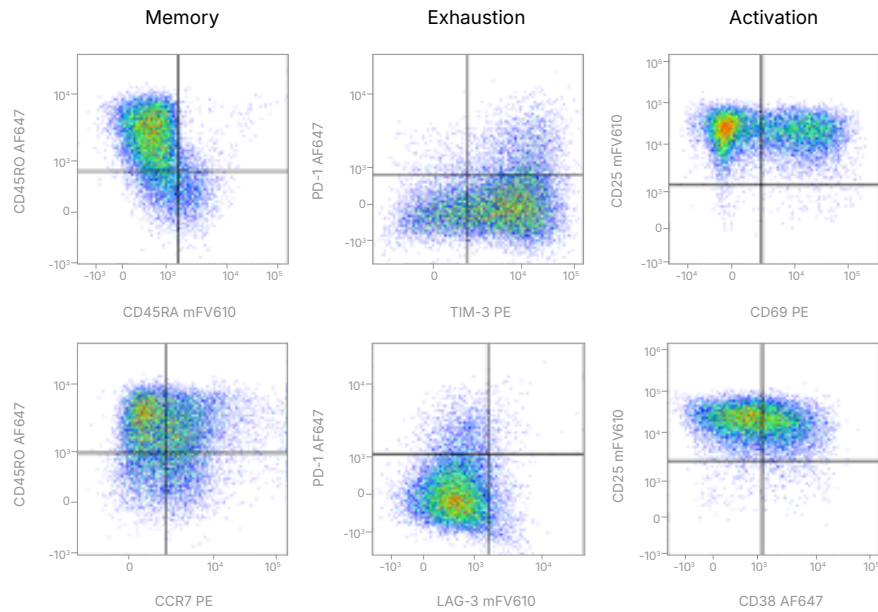


Figure 3: Characterization of Activated T Cells for T Cell Memory, Activation and Exhaustion with Multi-Color T-Cell Panels using Flow Cytometry. Day 2 activation and Day 9 memory and exhaustion phenotypes of expanded T cells. Human PBMCs stimulated with biotinylated versions of GMP Anti-Human CD3 (Cat# MAB11411-GMP) and GMP Anti-Human CD28 (Cat# MAB11412-GMP) coupled to M-270 Dynabeads + 200 IU/mL GMP Recombinant Human IL-2 (Cat# BT-202-GMP) for 9 days. Day 9 expanded T cells were stained with Anti-Human CD3 mFluor™ Violet 450, CD4 mFluor™ Violet 500, CD8 Alexa Fluor® 700, and Live/Dead 750/777 viability dye, followed by staining with three different panels to assess T cell 1) Memory (CD45RO Alexa Fluor® 647, CD45RA mFluor™ Violet 610, and CCR7 PE), 2) Exhaustion (PD-1 Alexa Fluor® 647, TIM-3 PE, and LAG-3 mFluor™ Violet 610), and 3) Activation (CD25 mFluor™ Violet 610, CD69 PE, and CD38 Alexa Fluor® 647). Learn more about [GMP antibodies](#) for immune cell therapy.

Flow Cytometry Panels Can Detect Donor-to-Donor Variability in Memory Phenotype, Activation, and Exhaustion Markers

Figure 4 demonstrates how using flow panels from Day 0 to Day 2 (for activation) or Day 9 (for memory and exhaustion) can provide quantitative data on how T cells from donors can change over time.

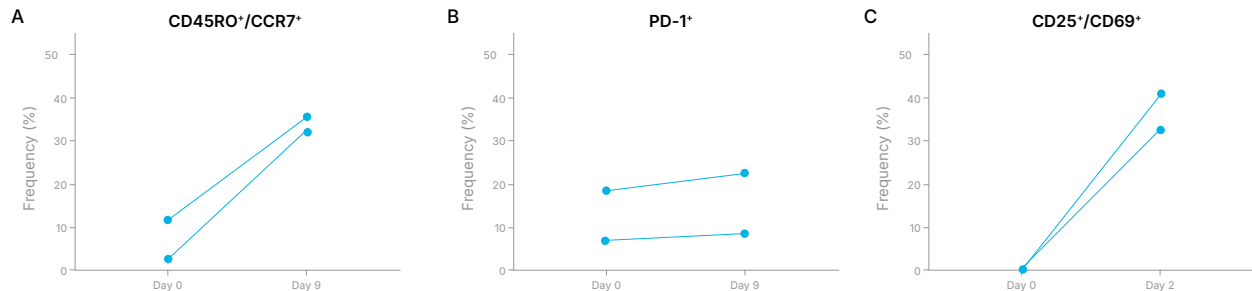


Figure 4: T Cell Memory, Activation and Exhaustion Panels Track Changes in T Cell Populations Over Time. Human PBMCs were stimulated with biotinylated versions of GMP Anti-Human CD3 (Cat# MAB11411-GMP) and GMP Anti-Human CD28 (Cat# MAB11412-GMP) coupled to M-270 Dynabeads + 200 IU/mL GMP Recombinant Human IL-2 (Cat# BT-002-GMP) for 9 days, as described in Figure 2. Differences in expression of CD45RO⁺/CCR7⁺ central memory (A), PD-1⁺ (B), and CD25⁺/CD69⁺ cells were analyzed at Day 0 and Day 9 (A, B), and Day 2 (C) after activation.

Multi-Color Antibody Panels are Compatible with Fluorokines for Flow Cytometry

Upon vector delivery and genetic engineering of the CAR-T cell, using flow panels to identify CAR-expressing T cells is necessary to ensure successful engineering of the cells and final product characterization before infusion back into the patient (Figure 1: Stage 4, Stage 6). Post-infusion of CAR-T cells, flow panels can be used to monitor disease progression, immune response, and the efficacy of CAR-T cell therapy (Figure 1: Stage 7).

Bio-Techne offers a wide selection of fluorescent-labeled recombinant proteins called Fluorokines™. Fluorokines are conjugated to Alexa Fluor® 488, Alexa Fluor® 647, Atto 488, or Atto 647N, offering intense fluorescence and excellent photostability. As demonstrated in Figure 5 below, Fluorokines can be combined with T cell antibody panels for comprehensive characterization of CAR-T cells using multi-color flow cytometry.

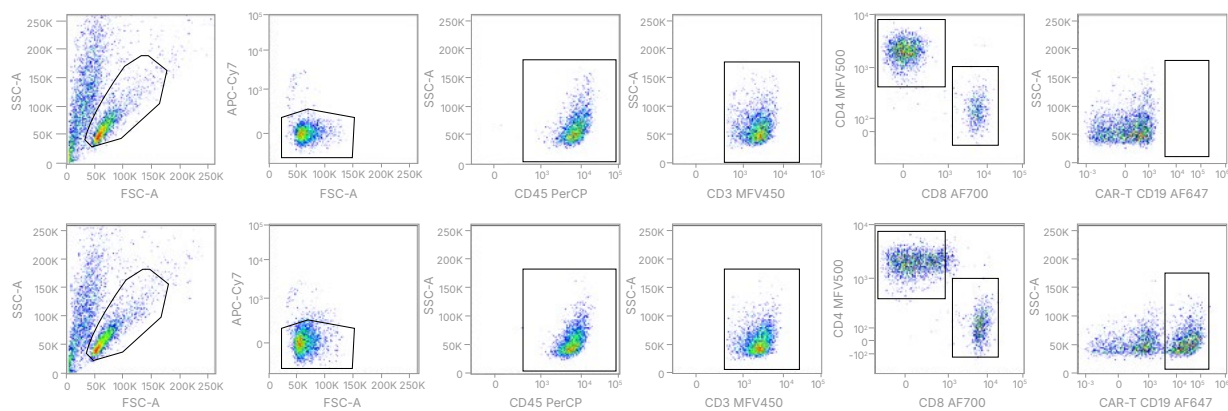


Figure 5. Characterization of CD4⁺/CD8⁺ CD19-CAR-T Cells by Multi-Color Flow Cytometry with a Recombinant Human CD19 CAR (AF647) Fluorokine and a Panel of Fluorochrome-Conjugated Antibodies. Two samples of CD4⁺/CD8⁺ T cells transduced with a hCD19-CAR and cultured were stained with the following panel of monoclonal antibodies: **PerCP-conjugated Mouse Anti-Human CD45** (R&D Systems, Cat# FAB1430C), **mFluor™ Violet 450-conjugated Mouse Anti-Human CD3 epsilon** (R&D Systems, Cat# FAB100MFV450), Alexa Fluor™ 700-conjugated **Mouse Anti-Human CD8 alpha** (R&D Systems, Cat# FAB1509N), **mFluor™ Violet 500-conjugated Mouse Anti-Human CD4** (R&D Systems, Cat# FAB3791MFV500), along with the Fluorokine, **Recombinant Human CD19 Fc Chimera Alexa Fluor® 647 Protein** (R&D Systems, Cat# AFR9269). Live/dead cells were identified using APC-Cy7. Cells were initially gated on singlets and live cells.

Conclusions

Bio-Techne's portfolio of immune cell marker antibodies enables multi-color T cell panels for comprehensive T Cell Memory, Activation, and Exhaustion immunophenotyping. These multi-color flow cytometry panels are compatible with our Fluorokines fluorescent-labeled proteins products, enabling detailed characterization of CAR-T cells. The flow cytometry panels used in this application note

show that Bio-Techne's R&D Systems antibodies can be used to immunophenotype naive PBMCs, activated PBMCs, and engineered CAR-T cells. All Bio-Techne reagents, including flow cytometry antibodies and Fluorokines, are backed by our 100% Guarantee. Bio-Techne ensures confidence in your results and the reliability of your data.

For research use or manufacturing purposes only. Trademarks and registered trademarks are the property of their respective owners. 6603962384

biotechne // Global Developer, Manufacturer, and Supplier of High-Quality Reagents, Analytical Instruments, and Precision Diagnostics.

INCLUDES R&D Systems™ Novus Biologicals™ Tocris Bioscience™ ProteinSimple™ ACD™ ExosomeDx™ Asuragen™ Lunaphore™