

# Exploring full-length T cell receptor sequencing and CITE-seq to identify and characterize pathogenic T cells in psoriasis

Ali Rezvan<sup>1</sup>, Cynthia Sakosfky<sup>1</sup>, Gracia Genero<sup>1</sup>, Lorena Hidalgo<sup>2</sup>, Hye-Won<sup>2</sup>, Aruna Ayer<sup>1</sup>, Stephanie Widmann<sup>2</sup>, Aaron J. Tzysnik, Gisele V. Baracho<sup>2</sup>

<sup>1</sup>BD Biosciences, Milpitas, California, USA <sup>2</sup>BD Biosciences, San Diego, California, USA,

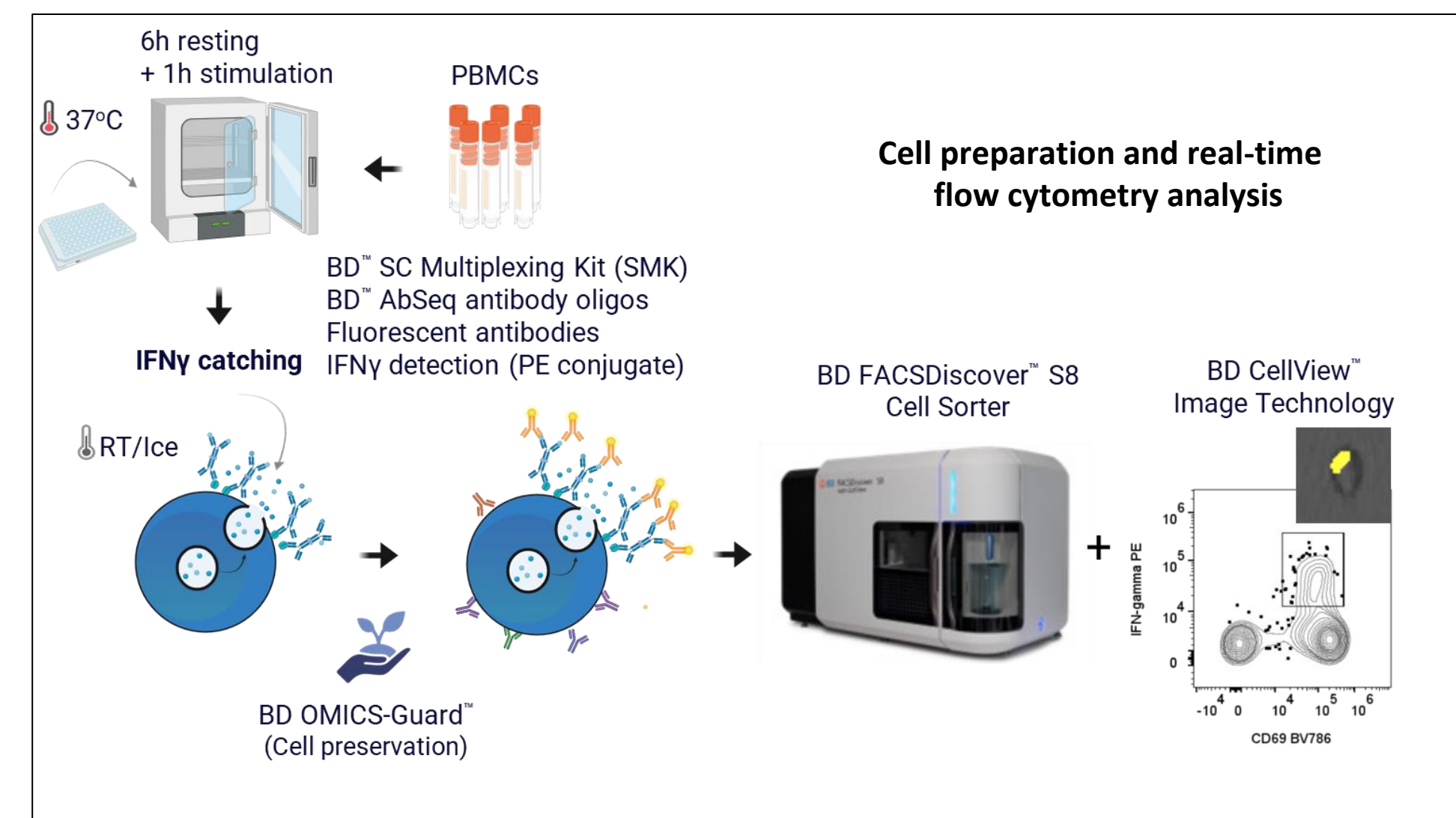
## Abstract

Psoriasis is a chronic immune-mediated skin disease characterized by keratinocyte hyperproliferation and systemic inflammation. In addition to its cutaneous manifestations, psoriasis is associated with comorbidities such as cardiovascular disease and psoriatic arthritis. The disease is driven by dysregulated T helper cells, which interact with dendritic cells and keratinocytes to sustain inflammation and recruit cytotoxic CD8<sup>+</sup> T cells. Early in disease progression, memory T cells, expressing cutaneous lymphocyte antigen (CLA) and skin-homing chemokine receptors, migrate into the skin and circulate between skin and blood, serving as peripheral biomarkers of disease activity. CLA<sup>+</sup> T cells, while less involved in lesion formation, may contribute to disseminating inflammation beyond the skin. To investigate the relationship between CLA<sup>+</sup> and CLA<sup>-</sup> T cells, we applied a multiomic approach combining single-cell full-length TCR sequencing, CITE-seq, and imaging flow cytometry. Imaging flow cytometry enabled real-time assessment of IFN $\gamma$  secretion without compromising cell viability. TCR sequencing revealed 263 shared clonotypes between CLA<sup>+</sup> and CLA<sup>-</sup> CD8<sup>+</sup> memory T cells, suggesting common antigen recognition. CLA<sup>+</sup> T cells from psoriatic donors were hyporesponsive and exhibited transcriptional exhaustion. In contrast, CLA<sup>-</sup> CD8<sup>+</sup> T cells secreted IFN $\gamma$  upon stimulation and showed a transcriptional program marked by expression of type I IFN genes and capable of rescuing these cells from exhaustion. This phenotype differed from healthy controls, where CLA<sup>-</sup> CD8<sup>+</sup> T cells displayed canonical cytotoxic profiles. These findings highlight the complementary roles of CLA<sup>+</sup> and CLA<sup>-</sup> T cells in local and systemic inflammation and demonstrate the utility of integrated multiomic technologies for understanding immune cell dynamics and accelerating biomarker discoveries in psoriasis.

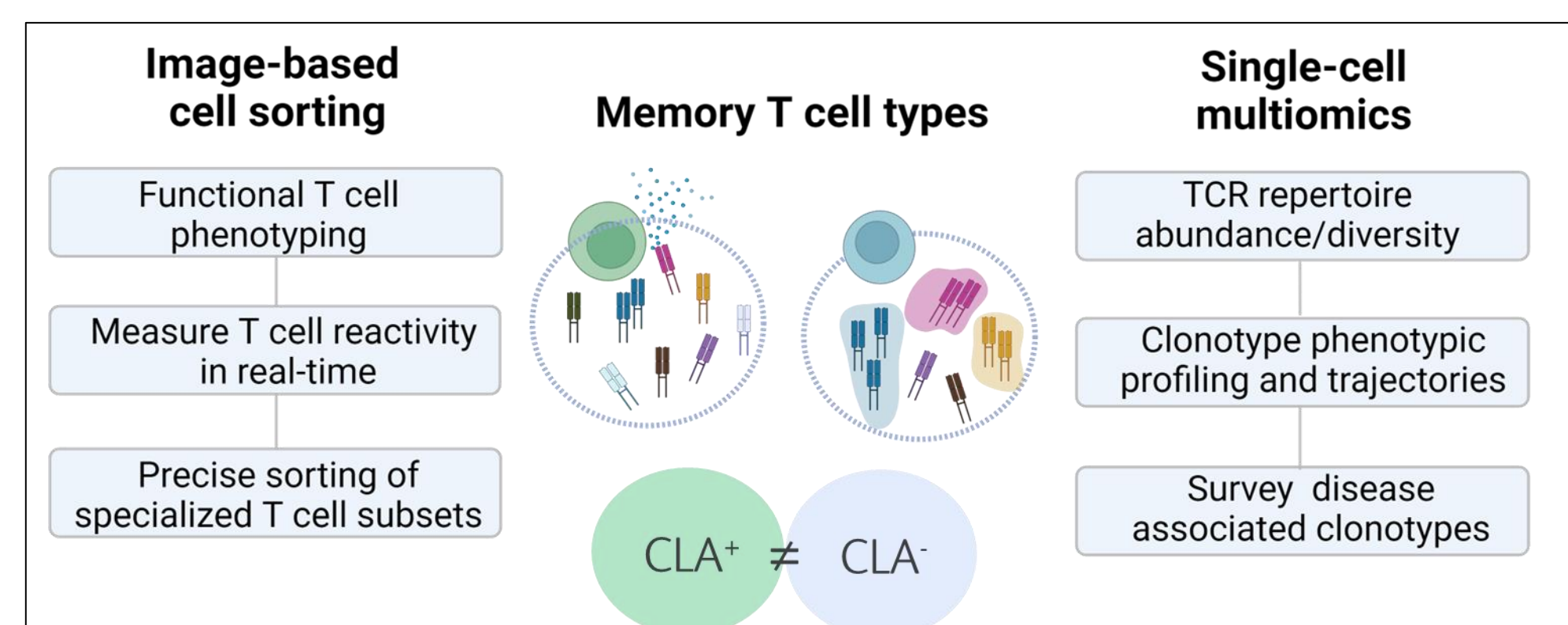
## Methods

### Integration of image-based cell sorting and single-cell multiomics for assessment of T-cell driven inflammation in psoriasis

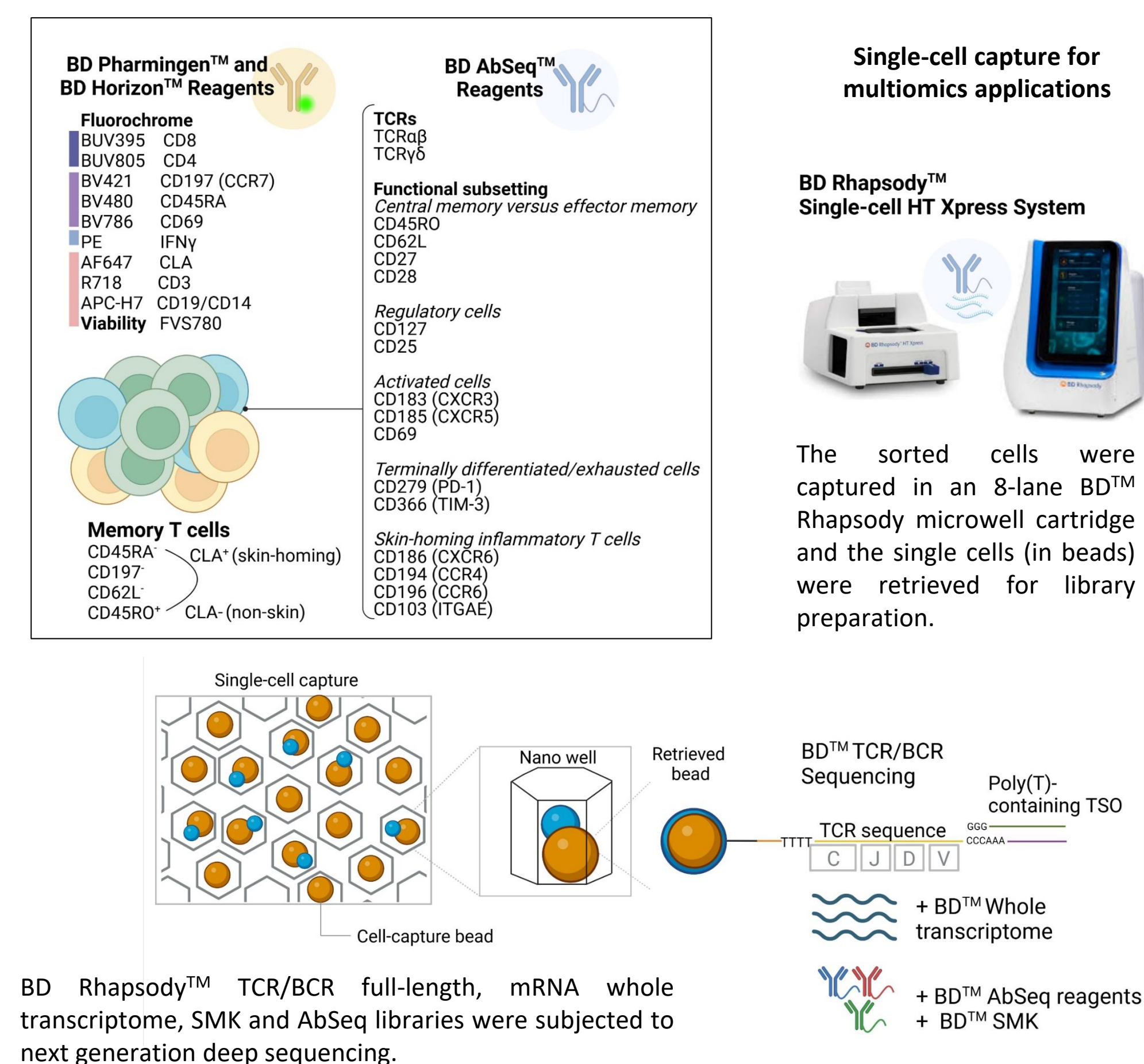
Frozen peripheral blood mononuclear cells from three healthy and three psoriatic individuals were incubated for 1 hour with CytoStim (Myltenyi Biotech) for restimulation of effector/memory T cells. After stimulation, secreted IFN $\gamma$  was measured using the Human IFN $\gamma$  Secretion Assay (Myltenyi Biotech) and the BD FACSDiscov<sup>TM</sup> S8 Imaging Cell Sorter.



In addition, cells from the different donors and conditions were barcoded with BD<sup>®</sup> Single Cell Multiplexing Kit tags and simultaneously labeled with antibody-oligos and fluorescent antibodies. This enabled the sorting of memory subsets for downstream single-cell TCR sequencing and CITE-seq analyses.

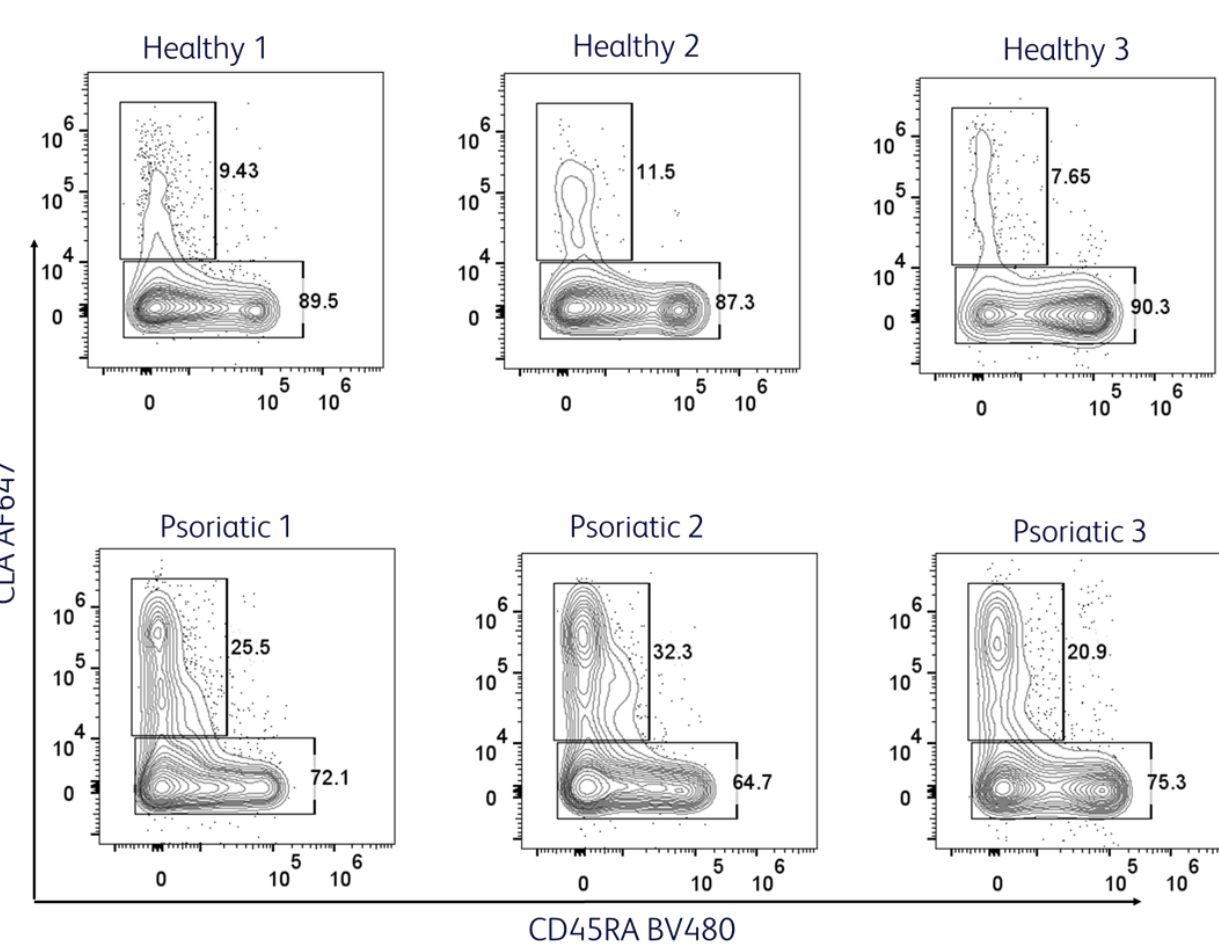


Antibody panels were designed for characterization of memory T cell subsets, including skin-derived cutaneous lymphocyte antigen (CLA)<sup>+</sup> T cells. The sorted populations CD3<sup>+</sup>CD45RA<sup>+</sup>CD197<sup>+</sup>CLA<sup>+</sup> or CD3<sup>+</sup>CD45RA<sup>+</sup>CD197<sup>+</sup>CLA<sup>-</sup> were obtained from both unstimulated and stimulated cell cultures.

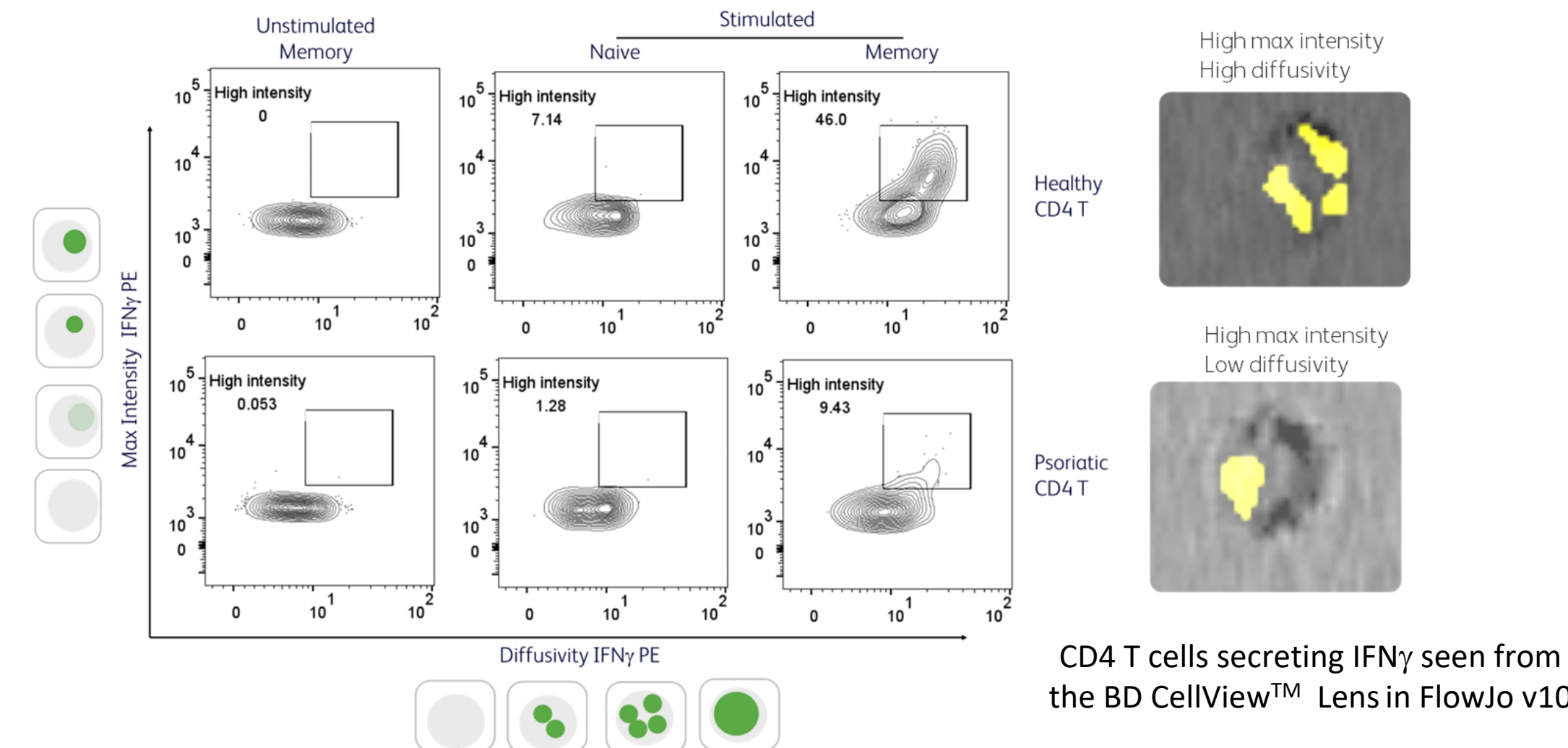


## Results – Functional fitness of circulating skin-homing T cells

### 1A High frequencies of skin-homing CLA<sup>+</sup> CD4T cells in psoriasis



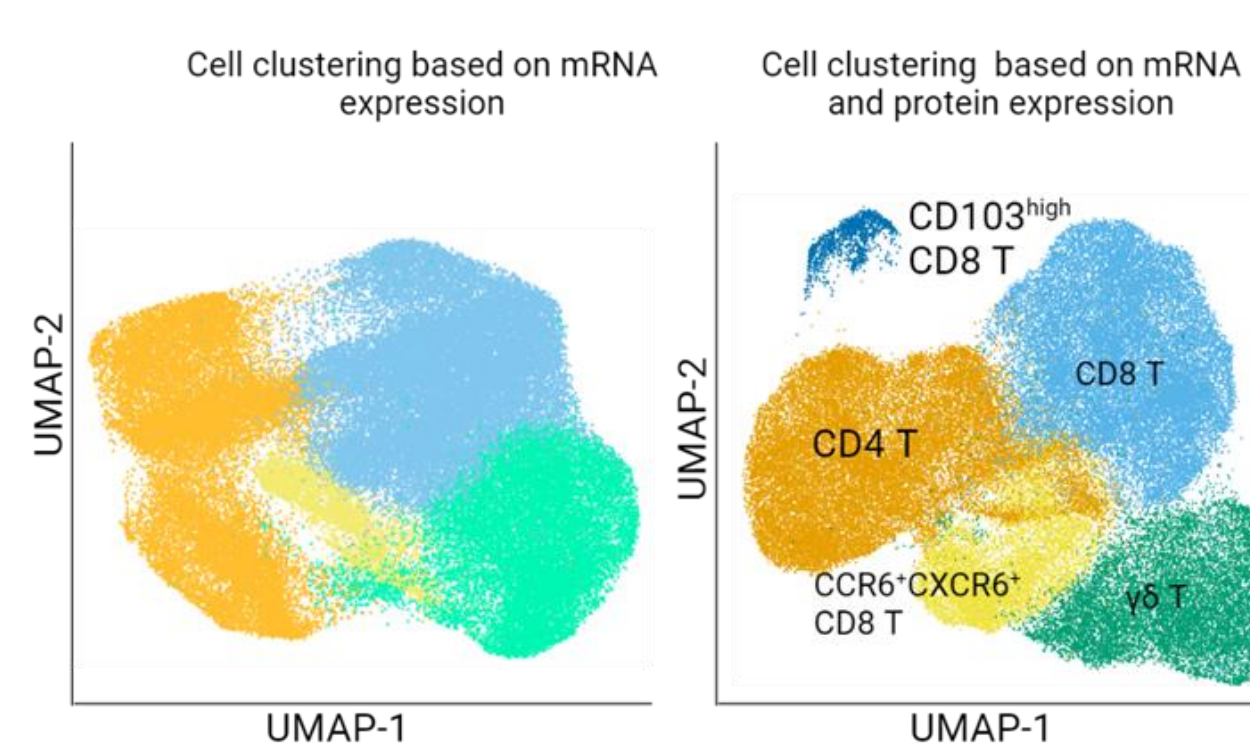
### 1B Defective IFN $\gamma$ secretion by CD4T cells but not CD8 T cells in psoriasis



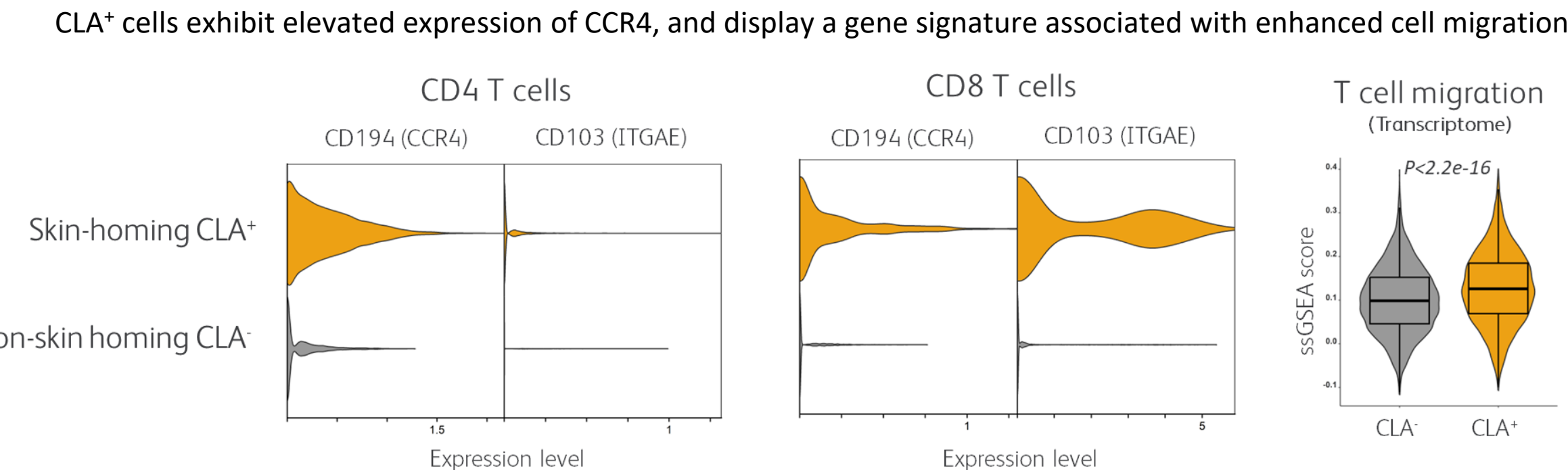
**Figure 1: Functional fitness of circulating T cells measured by imaging flow cytometry.** A) Assessment of the frequencies of CLA<sup>+</sup> T cells among PBMCs. B and C) Detection of functional T cells secreting IFN $\gamma$  using imaging fluorescence parameters on the BD FACSDiscov<sup>TM</sup> S8 Cell Sorter upon re-stimulation of memory T cells with CytoStim. C) Further assessment of IFN $\gamma$  secretion reveals that CLA<sup>+</sup> cells are less responsive to CytoStim stimulation and unlikely secrete IFN $\gamma$ .

## Results – Proteomic and transcriptomic characteristics of skin-homing T cells

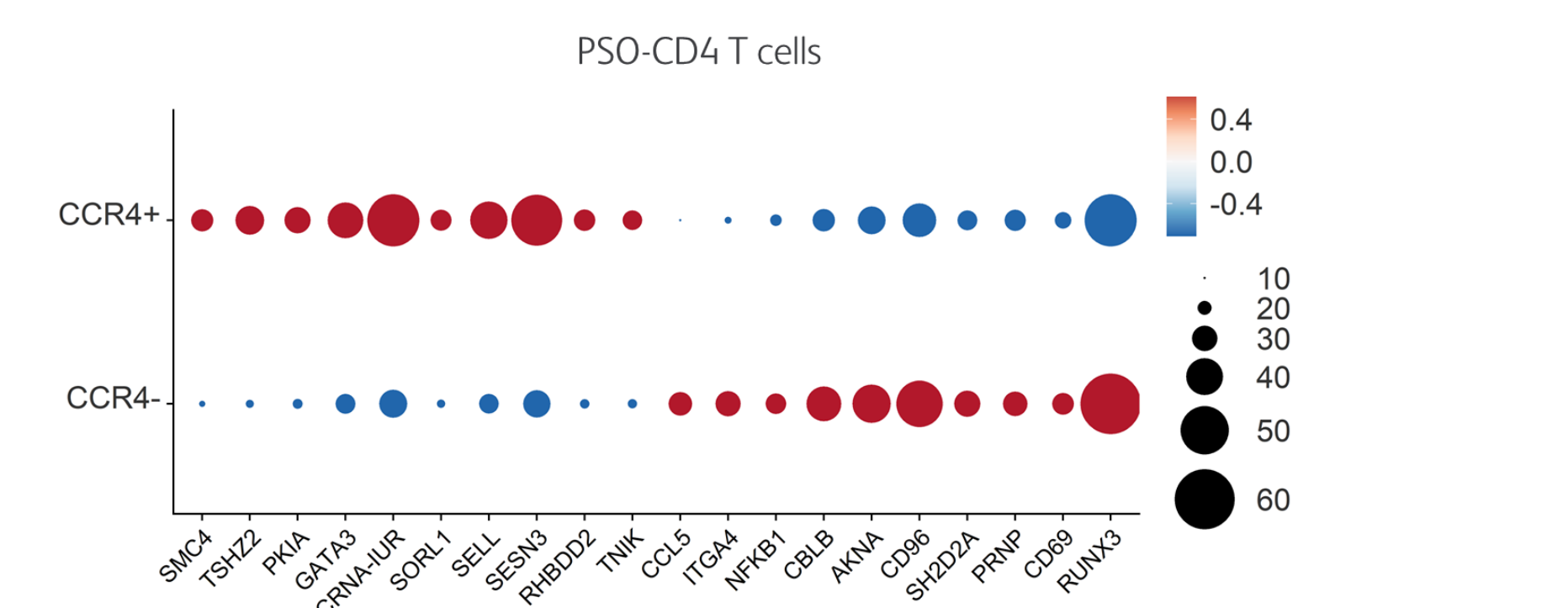
### 2A Optimized detection of circulating CD103<sup>+</sup> cells with CITE-seq



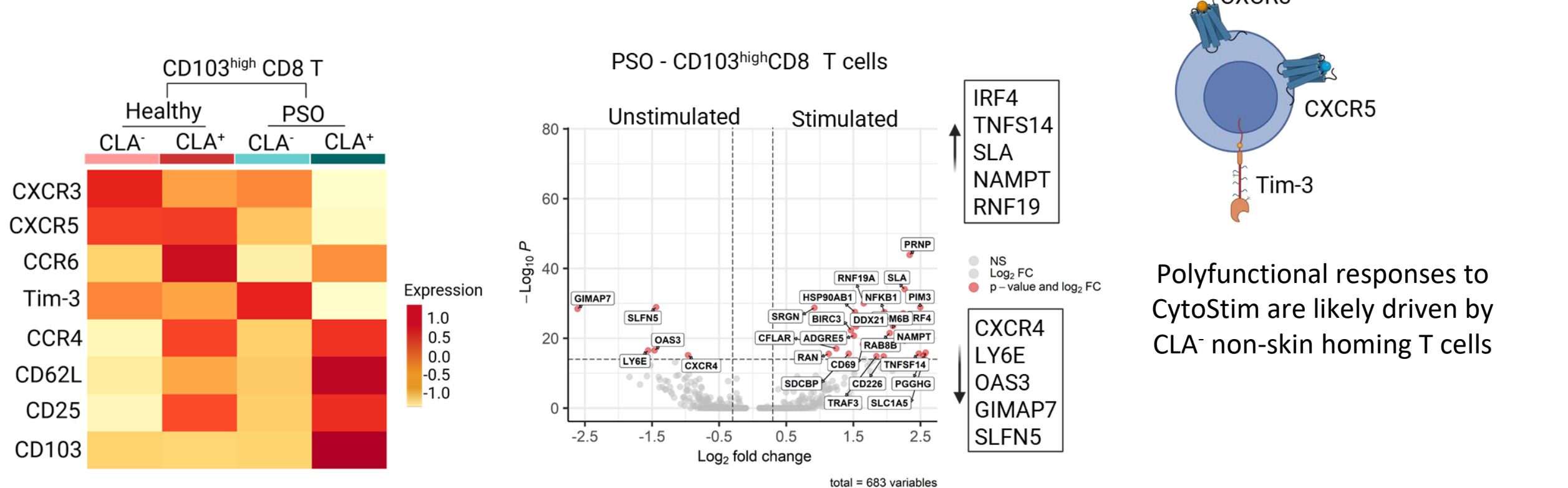
### 2B Skin-homing T cells are enriched for cell migration-related genes



### 2C Skin-homing CCR4<sup>+</sup> CD4 T cells from psoriatic individuals exhibit a gene expression profile consistent with a senescent/exhausted phenotype

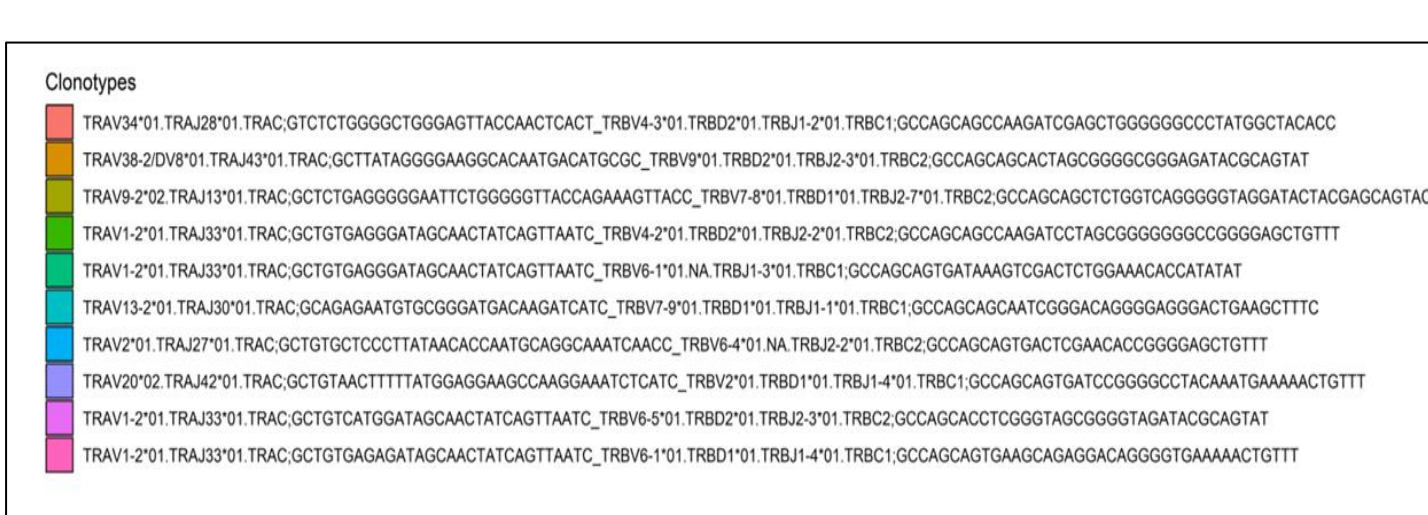
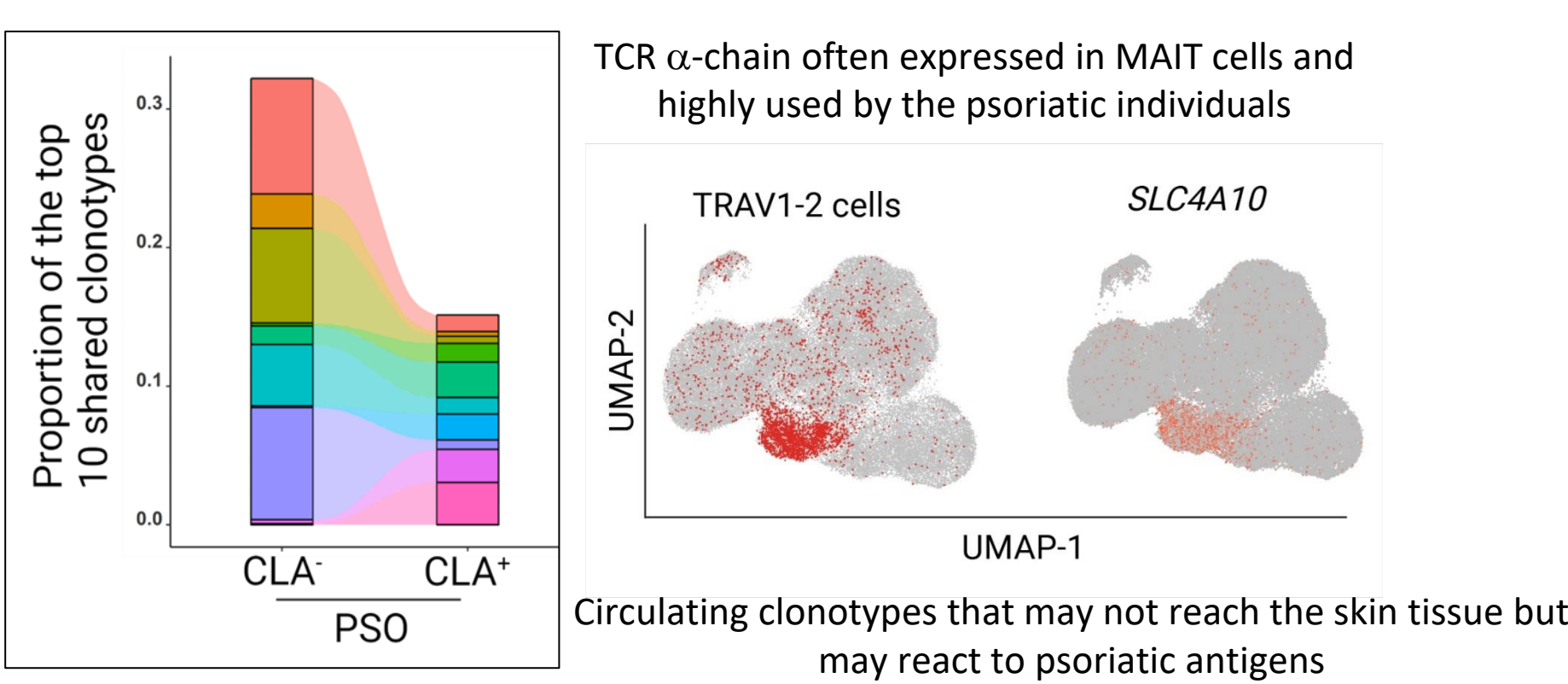


### 2D CD103<sup>+</sup> CD8 T cells from psoriatic individuals exhibit impaired expression of key effector markers



**Figure 2: Phenotypic and functional profiling of skin-homing and non-skin homing memory T cells using CITE-seq.** A) Dimensionality reduction and Seurat clustering of over 60,000 high quality cells demonstrate improved spatial cell organization when clustering is performed using integrated mRNA and protein data. B) Gene enrichment analysis reveals that skin-homing cells exhibit higher score for cell migration gene signatures compared to non-skin homing cells, supporting their ability to migrate between skin tissue and blood. C) Differential gene expression analysis of CCR4<sup>+</sup> and CCR4<sup>-</sup> CD4 T cells from psoriatic individuals identifies elevated expression of *SESN3*, *LCNRA-IUR* and *TSHZ2* that may contribute to T cell exhaustion and impaired IFN $\gamma$  secretion in response to CytoStim. D) Protein analysis highlights defective expression of CXCR3, CXCR5 and Tim-3 in CLA<sup>+</sup> skin-homing T cells from psoriatic individuals. Despite lower expression of these effector targets compared to healthy individuals, the CD103<sup>+</sup>CD8 T cells from the psoriatic individuals retain responsiveness to CytoStim stimulation. This response is mostly driven by the CLA<sup>+</sup> cells counterpart that express effector protein markers

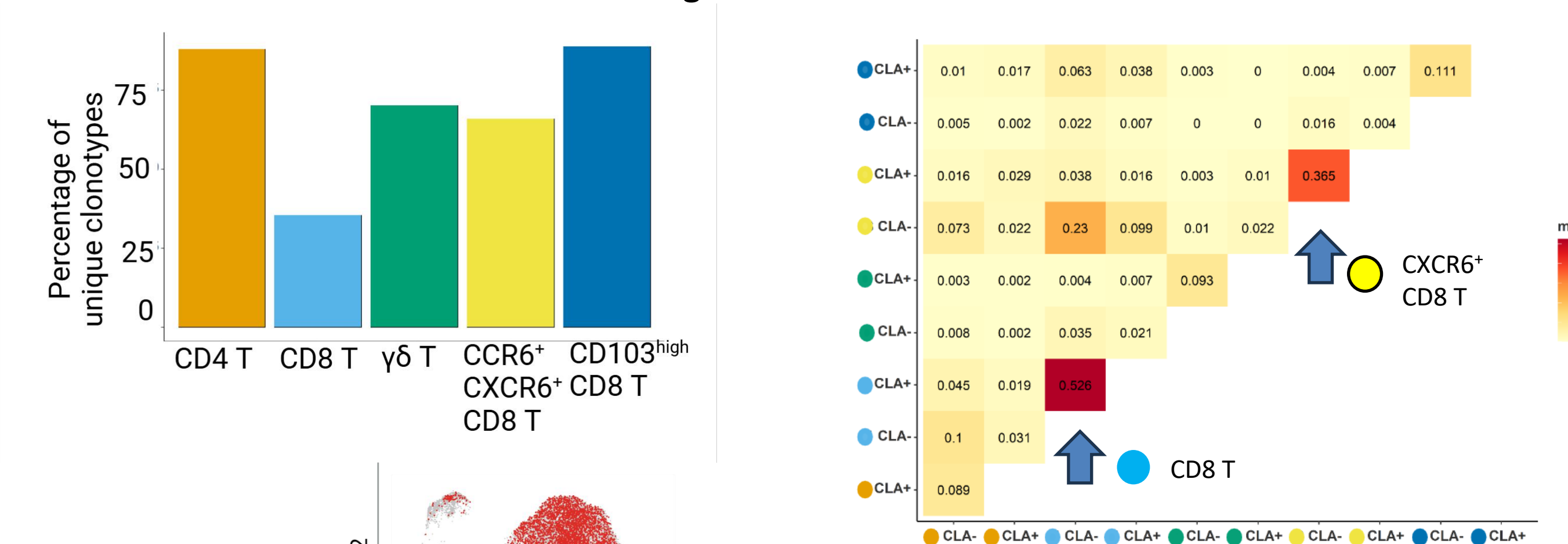
### 3B Skin- and non-skin-homing CD8 T cells can share TCR clonotypes while exhibiting distinct functional profiles



**B) The alluvial plot depicts the top 10 clonotypes shared between skin- and non-skin homing T cells, each identified by a distinct color. The UMAP displays one of the most frequently used  $\alpha$ -chain genes in the psoriatic individuals, which is also shared with non-skin homing cells. The shared clones from healthy individuals in addition to having similar TCR sequences, exhibit overall similar functional phenotype. In contrast, the shared clones from psoriatic individuals present distinct transcriptional programs: CLA<sup>+</sup> cells are suppressed and become exhausted upon stimulation, while CLA<sup>-</sup> cells retain INF I responses.**

## Results – Full-length TCR sequencing uncovers clones shared between skin- and non-skin homing T cells

### 3A Non-skin homing CD8 T cell subsets share TCR clonotypes with skin-homing CD8 T cells



## Conclusions

Using imaging flow cytometry, we assessed the functional fitness of skin-homing T cells through real-time analysis of IFN $\gamma$  secretion. Additionally, we sorted memory subsets for comprehensive downstream multiomics analysis using a complete workflow solution provided by the BD<sup>®</sup> Rhapsody Single Cell Capture System.

This workflow eliminated the need for skin tissue preparations in analyzing skin-derived T cells and further enabled the identification of TCR clonotypes shared between peripheral memory and skin-homing memory T cells.

In the psoriatic samples, CLA<sup>+</sup> cells expressing overlapping clonotypes with skin-homing cells upregulated *IRF4*, *PRDM1* and other genes suggesting that these cells can respond to stimulation and sustain type I IFN-mediated inflammation during disease flares. Importantly, these cells may encounter specific antigens outside the skin, contributing to spread of the disease to other locations.

In contrast, CLA<sup>-</sup> skin-homing cells exhibited signs of suppression and may undergo exhaustion upon stimulation.

Our results help to elucidate immune mechanisms involved in psoriasis, offering insights that may inform future research into targeted treatments for this and related skin inflammatory conditions.

BD Flow Cytometers are Class I Laser Products. For Research Use Only. Not for use in diagnostic or therapeutic procedures. BD and the BD logo are trademarks of Becton, Dickinson and Company. All other trademarks are the property of their respective owners. © 2025 BD. All rights reserved. BD-161583 (v1.0) 0925