

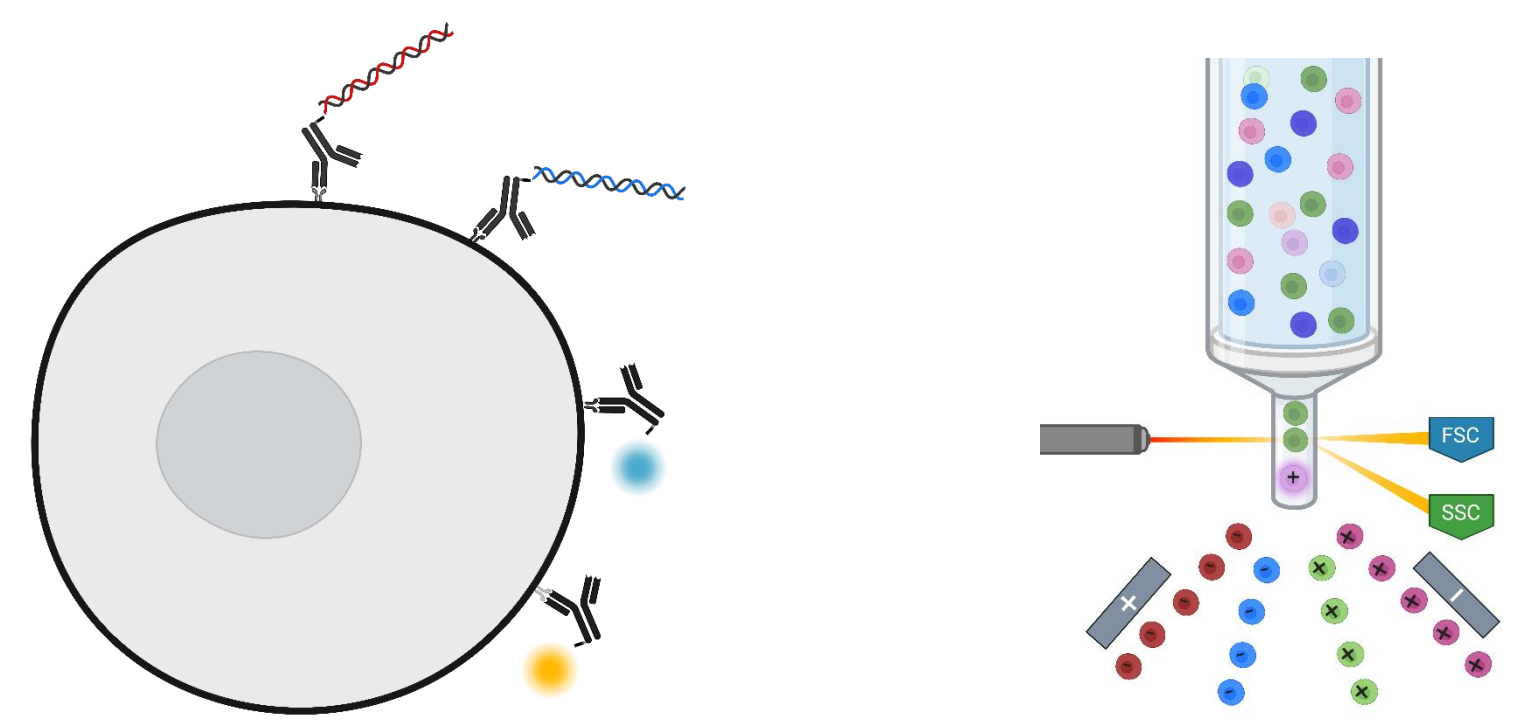
Abstract

T follicular helper (T_{FH}) cells are a CD4⁺ T cell subset that play a critical role in the formation of germinal centers (GC). T_{FH} cells are essential to produce protective antibody responses through the regulation of clonal selection and differentiation into activated antibody-secreting and memory B cells. T_{FH} cell differentiation is a multistep process that begins with the presentation of antigen by dendritic cells to naïve CD4⁺ T cells, where the CD4⁺ T cells commit to the T_{FH} lineage by expressing the lineage defining transcription factor B cell lymphoma 6 (Bcl-6). T_{FH} cells are phenotypically characterized by the high expression of C-X-C chemokine receptor 5 (CXCR5) and programmed cell death-1 (PD-1) on the cell surface. CXCR5⁺ memory CD4⁺ T cells present in peripheral blood represent a circulating T_{FH} (cT_{FH}) cell subset. In contrast to T_{FH} cells, cT_{FH} cells may express low levels of Bcl-6 with numerous reports of subsets that differ both phenotypically and functionally.

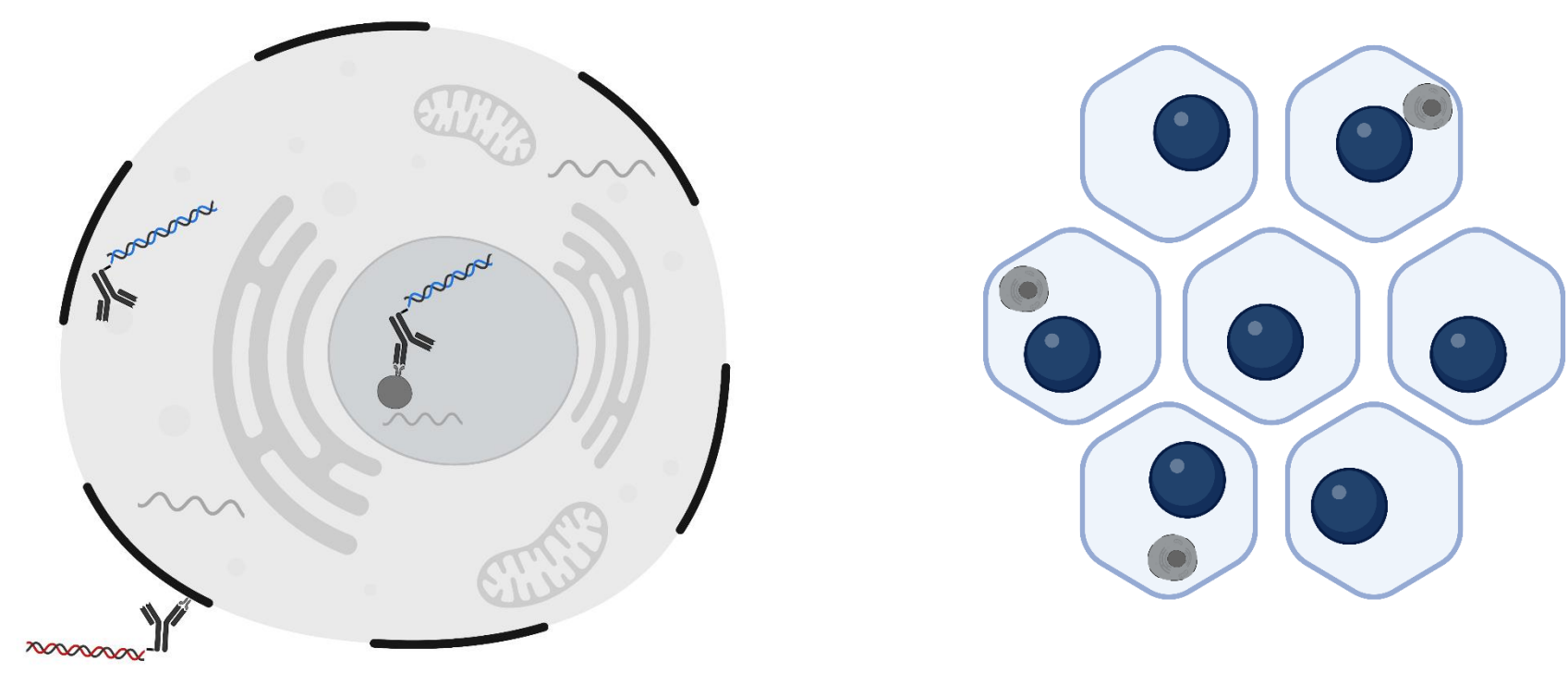
Here we characterize the heterogeneous subsets of T_{FH} and cT_{FH} cells using a combination of intracellular CITE-seq, surface CITE-seq and whole transcriptome single-cell RNA sequencing (scRNA-seq) analysis. Germinal center and circulating T_{FH} cells were enriched by sorting memory CD4⁺, CXCR5⁺, PD-1⁺ from a tonsil (n=1), and memory CD4⁺ CXCR5⁺ cells from a matched blood sample and a healthy blood donor (n=2). We examined the expression of Bcl-6 and other transcription factors, using protein and RNA expression data, across pre-T_{FH} (an early development stage in the germinal center), T_{FH} and cT_{FH} cells. Germinal center and circulating T_{FH} cells profiles were also compared.

Together, these data demonstrate how a combination of intracellular CITE-seq, surface CITE-seq and scRNA-seq may be used to deepen our understanding of the differentiation states and subsets of cells present at low abundance, such as T_{FH} cells, by enabling detection of intracellular and surface protein along with high-quality mRNA.

BD Rhapsody™ System Workflow



Single-cell suspensions from tonsil and blood were stained with a panel of antibodies including CD4, CD45-RA, PD-1 and CXCR5 to identify T_{FH} cells. Cells were also labeled with a 23-plex surface BD[®] AbSeq Panel. Cells were sorted using a 4-way sort on the BD FACSMelody™ Cell Sorter.



T_{FH} enriched immune cell suspensions were fixed using BD[®] Sample Preservation Buffer followed by permeabilization with BD Phosflow™ Perm Buffer III and stained with intracellular, oligo-conjugated BD[®] AbSeq Antibodies. Cells were counted with the BD Rhapsody™ System and loaded onto the cell capture cartridge. Libraries were prepared and indexed for Illumina sequencing followed by processing using the BD Rhapsody™ Whole Transcriptome Analysis Pipeline on Seven Bridges. Read count matrices were analyzed and visualized using Seurat and scCustomize in RStudio.^{1,2} Briefly, standard QC metrics were applied followed by dataset integration, normalization, data scaling and visualization.

T_{FH} Enrichment

Figure 1. The BD FACSMelody™ Cell Sorter sorting strategy

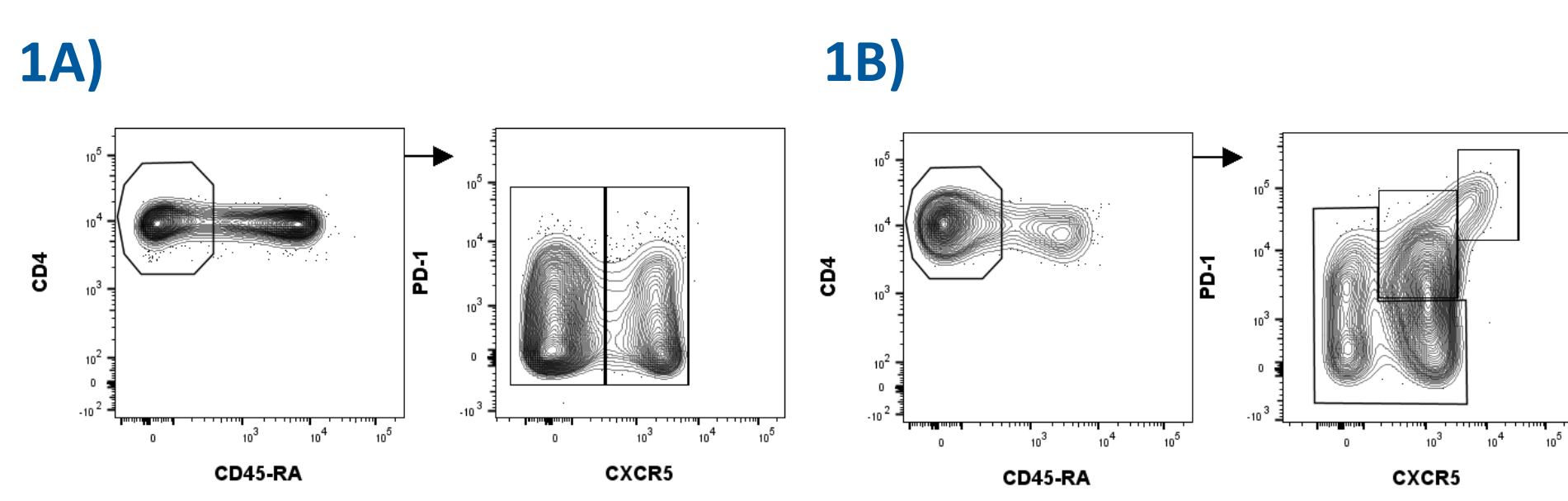


Figure 1. The BD FACSMelody™ Cell Sorter sorting strategy for 1A) cT_{FH} in blood and 1B) T_{FH} in tonsil prior to BD Rhapsody™ whole transcriptome workflow.

Identification and characterization of T_{FH} and cT_{FH} using intracellular and surface BD[®] AbSeq Antibodies

Figure 2. Tonsillar T_{FH} Cells

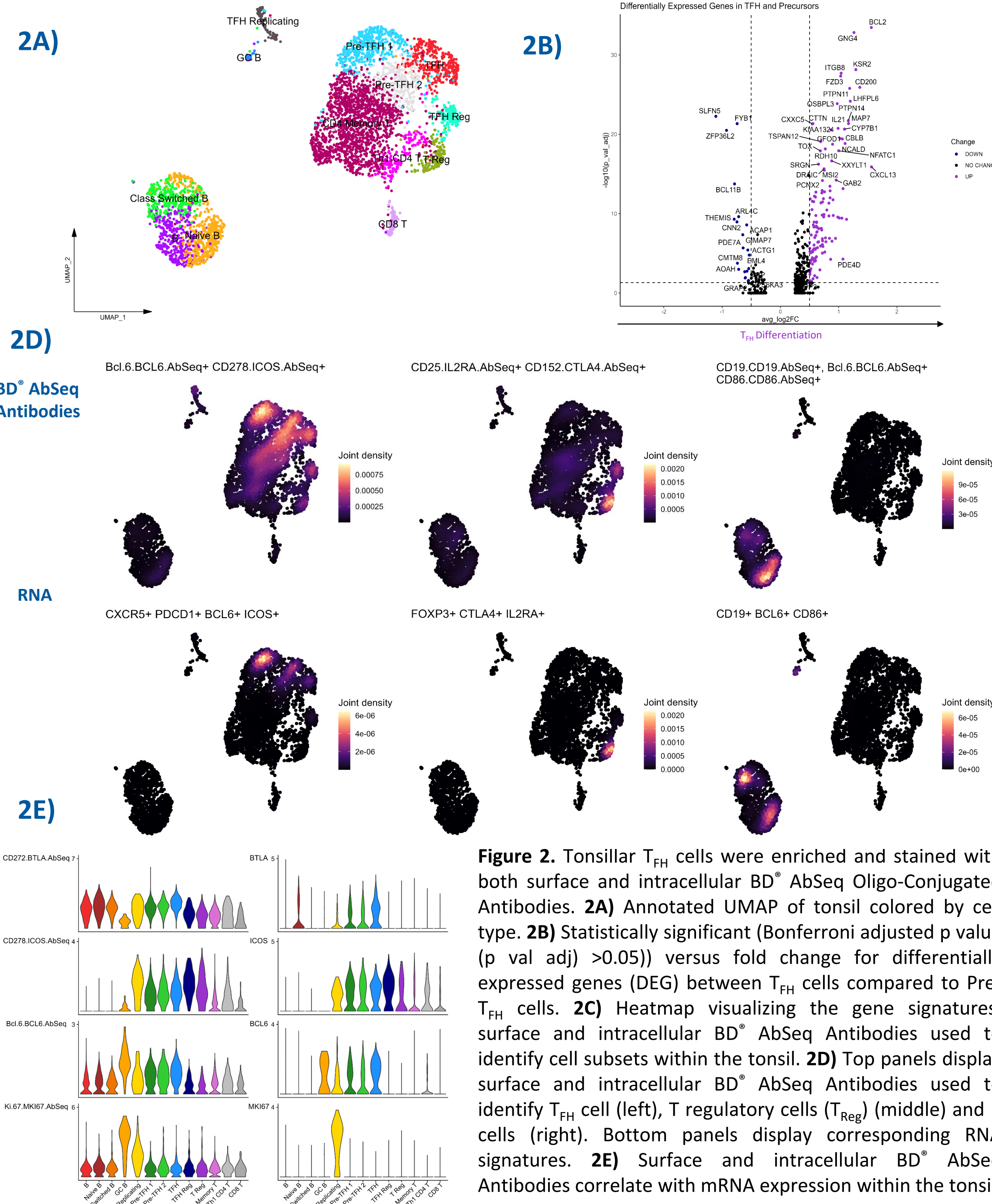


Figure 2. Tonsillar T_{FH} cells were enriched and stained with both surface and intracellular BD[®] AbSeq Oligo-Conjugated Antibodies. 2A) Annotated UMAP of tonsil colored by cell type. 2B) Statistically significant (Bonferroni adjusted p value (p val adj) >0.05) versus fold change for differentially expressed genes (DEG) between T_{FH} cells compared to Pre-T_{FH} cells. 2C) Heatmap visualizing the gene signatures, surface and intracellular BD[®] AbSeq Antibodies used to identify cell subsets within the tonsil. 2D) Top panels display surface and intracellular BD[®] AbSeq Antibodies used to identify T_{FH} cell (left), T regulatory cells (T_{Reg}) (middle) and B cells (right). Bottom panels display corresponding RNA signatures. 2E) Surface and intracellular BD[®] AbSeq Antibodies correlate with mRNA expression within the tonsil.

Figure 3. cT_{FH} Cells in Blood

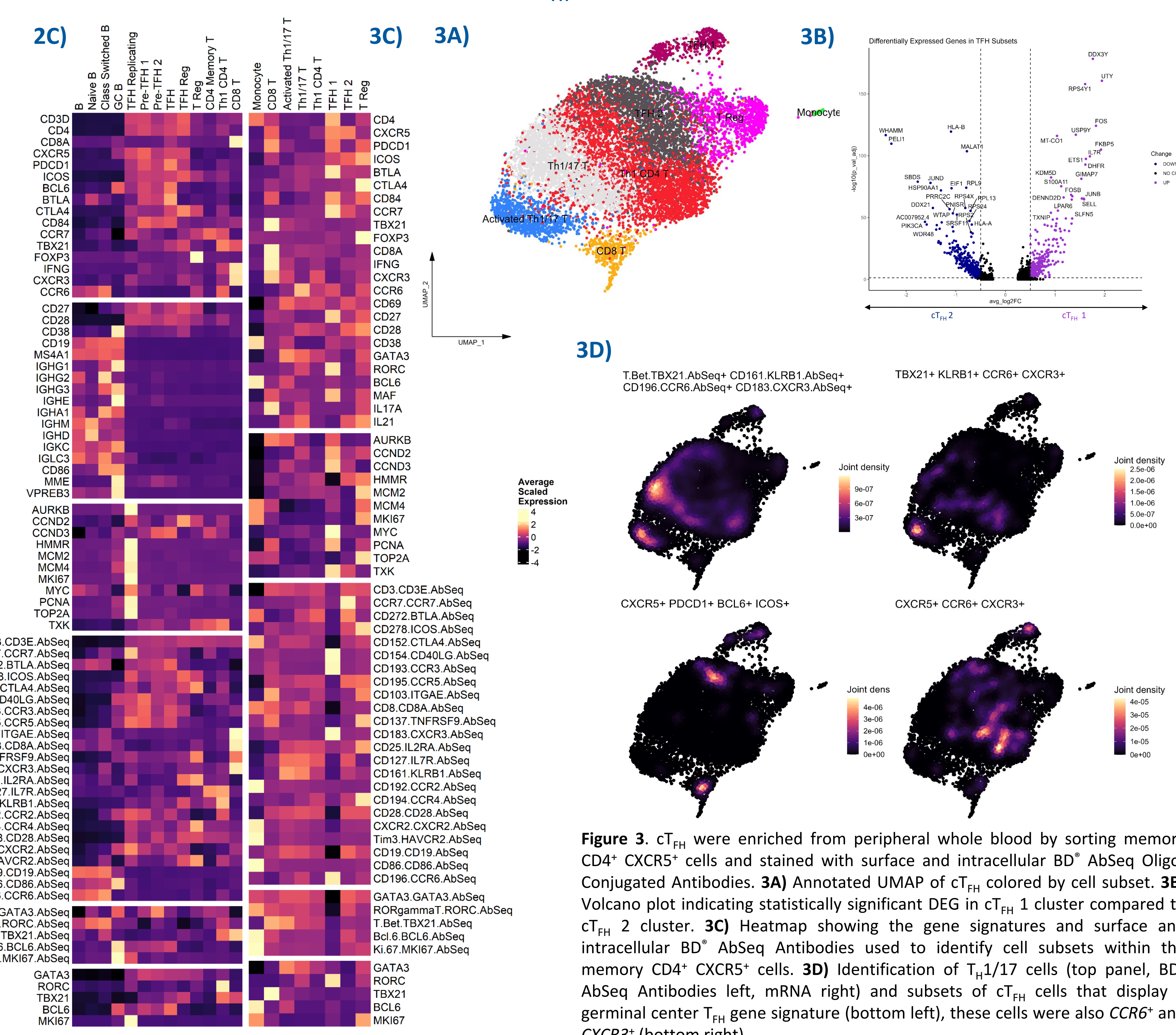


Figure 3. cT_{FH} were enriched from peripheral whole blood by sorting memory CD4⁺ CXCR5⁺ cells and stained with surface and intracellular BD[®] AbSeq Oligo-Conjugated Antibodies. 3A) Annotated UMAP of cT_{FH} cells colored by cell subset. 3B) Volcano plot indicating statistically significant DEG in cT_{FH} 1 cluster compared to cT_{FH} 2 cluster. 3C) Heatmap showing the gene signatures and surface and intracellular BD[®] AbSeq Antibodies used to identify cell subsets within the memory CD4⁺ CXCR5⁺ cells. 3D) Identification of T_{1H}/17 cells (top panel), BD[®] AbSeq Antibodies left, mRNA right) and subsets of cT_{FH} cells that display a germinal center T_{FH} gene signature (bottom left), these cells were also CCR6⁺ and CXCR3⁺ (bottom right)

Analysis of cell-cell communication

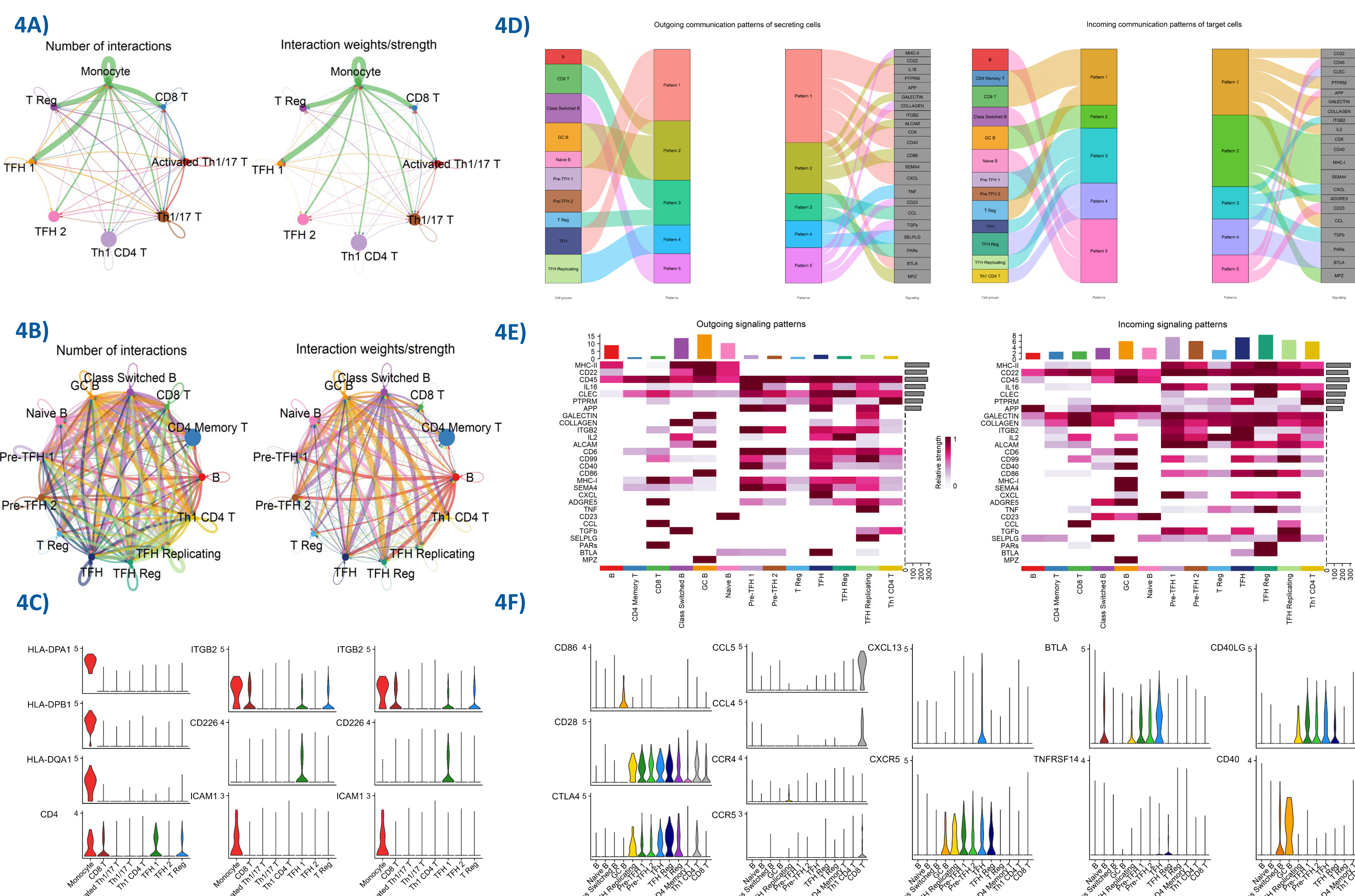


Figure 4. The probability of cell-cell communication was computed for tonsillar T_{FH} and cT_{FH} cells using CellChat.³ 4A) Aggregated cell-cell communication network with number of interaction and interaction weights for peripheral blood resident cells. 4B) Aggregated cell-cell communication network with number of interaction and interaction weights for tonsil resident cells. 4C) Gene expression of selected enriched signaling pathways in peripheral blood. 4D) The top outgoing and incoming signaling patterns of secreting cells and target cells visualized by river plot, which indicates the coordination of cell types to global tonsil function. 4E) Signals that contribute most to outgoing and incoming signals of cell groups in tonsil visualized by heatmap. 4F) Gene expression of selected enriched signaling pathways in tonsil.

Comparison of T_{FH} and cT_{FH}

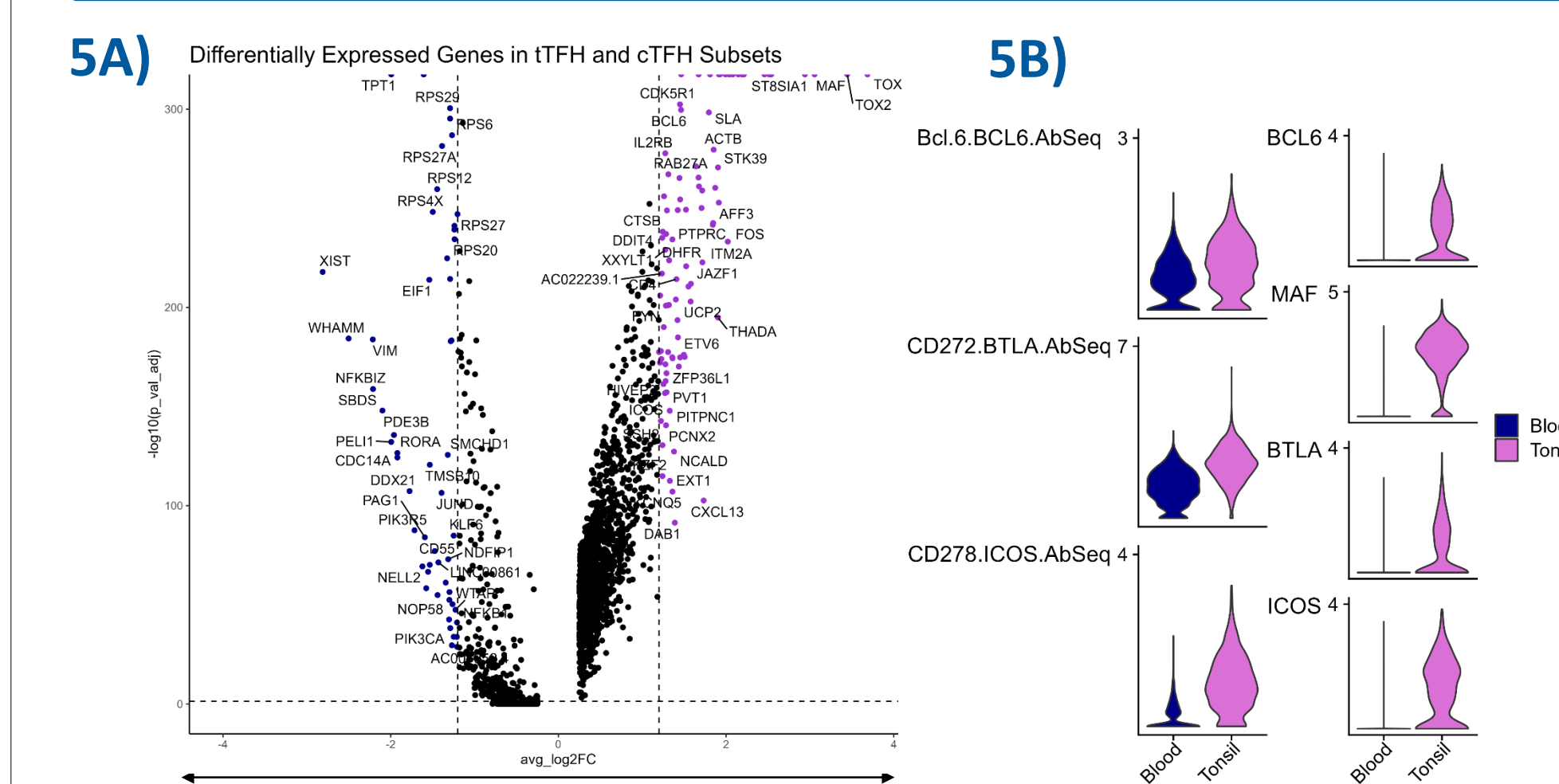


Figure 5. DEG were computed for tonsillar T_{FH} cells and cT_{FH}. 5A) Volcano plot indicating statistically significant DEG in tonsillar T_{FH} cells and cT_{FH}. 5B) Key gene and corresponding BD[®] AbSeq Antibody expression visualized with violin plot.

Conclusions

- We have demonstrated a workflow that enables the simultaneous detection of intracellular BD[®] AbSeq and surface BD[®] AbSeq Antibodies along with high-quality mRNA.
- Transcription factors detected with intracellular BD[®] AbSeq Antibodies showed a high concordance with mRNA expression.
- Our data demonstrate how a combination of intracellular BD[®] AbSeq Antibodies, surface BD[®] AbSeq Antibodies and scRNA-seq may be used to identify subsets of cells present at low abundance, such as T_{FH}, pre-T_{FH} and cT_{FH} cells.

References:
 1) Hao Y, Hao S, Andersen-Nissen E, et al. *Cell*. 2021;184(13):3573-3587.e29. doi: 10.1016/j.cell.2021.04.048.
 2) Marsh SE. 2021. doi: 10.5281/zenodo.5706430.
 3) Jin S, Guerrero-Juarez CF, Zhang L, et al. *Nat Commun*. 2021;12(1):1088. doi: 10.1038/s41467-021-21246-9.
 4) Some images were created with BioRender.com