

A fast and efficient bioinformatics analysis workflow for processing reads from single-cell multiomics assays captured on a microwell-based platform

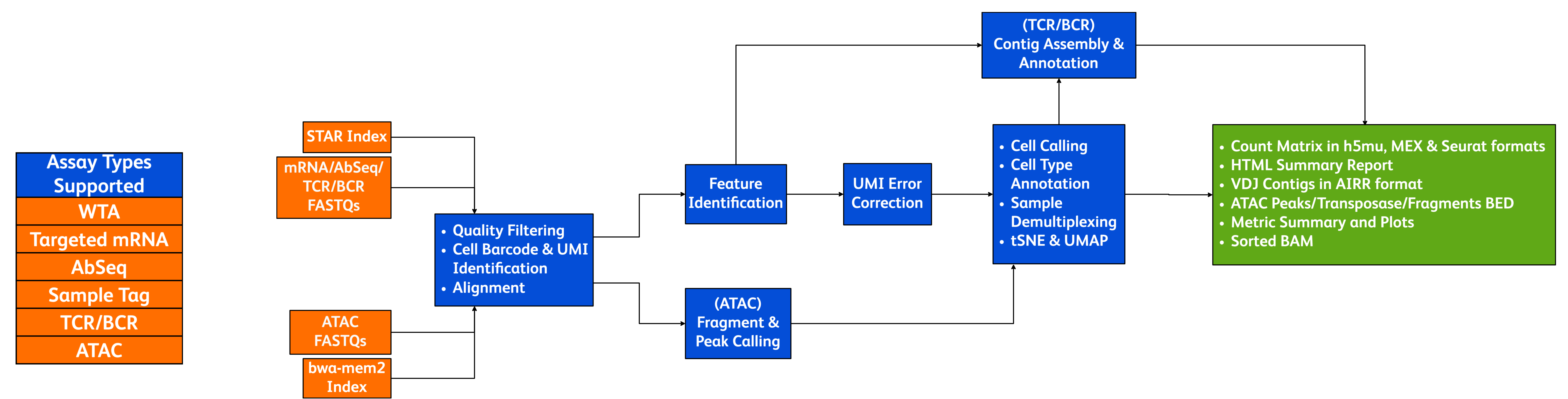
Poster# 4956

Raghavendra Padmanabhan¹, Brent Weichel¹, Amie Radenbaugh¹, Yuefu Jiang¹, Charles Weeks¹, Thomas McCarthy¹, Youngsook Kim¹, Anthony Berno¹, Devon Jensen¹.
¹BD Biosciences, 2350 Qume Drive, San Jose, CA 95131

Abstract

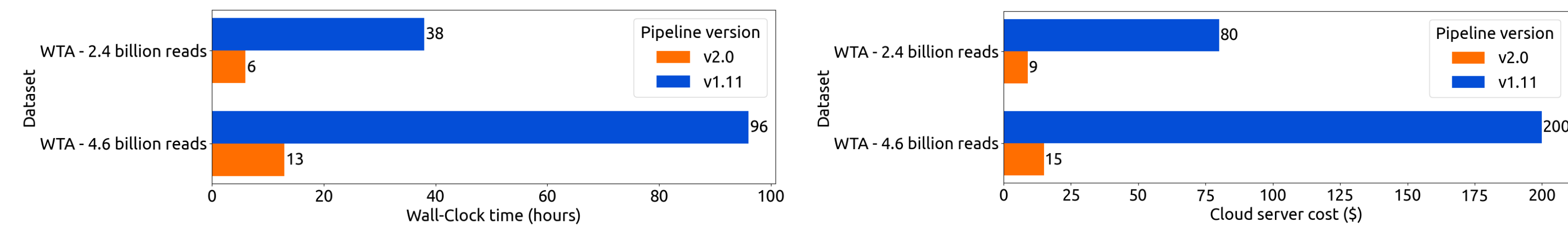
Here, we present a significant update of the BD Rhapsody™ Sequence Analysis Pipeline, version 2—a comprehensive primary analysis workflow that is up to 7X faster and consumes 2X less disk space than the previous release. We used it to process a library with 20,000 human PBMCs sequenced to a depth of 1.2 billion reads in 2.5 hours of wall-clock time. The pipeline can be used with reads from single-cell whole transcriptome, targeted mRNA, surface antigens (AbSeq), TCR/BCR, ATAC-seq and Sample Tag libraries captured on the BD Rhapsody™ System platform. The major processing steps include Quality Filtering, Cell Barcode Identification, Read Alignment, Feature Assignment, Fragment & Peak calling, UMI Error Correction, Identification of Putative Cells, Sample De-Multiplexing, Dimensionality Reduction for Visualization and VDJ Contig Assembly & Annotation. Output files and metrics are available in easy-to-digest formats, including an html report with dynamic visualization. Integrated outputs in Seurat and Scanpy formats combine expression matrices and all cell annotation metadata (e.g., predicted cell type, sample assignment, TCR/BCR sequence and gene segments). These pre-generated files are ready to load into popular single-cell analysis tools. The pipeline uses CWL as the workflow manager and makes use of custom code written in C++ and Python along with various open-source packages for data processing. The pipeline can also process reads from non-human species and includes built-in support for specifying and building custom reference genome indices. In support of high-throughput multiomic discovery studies enabled by the BD Rhapsody™ HT Xpress System, this pipeline has been tested with datasets containing more than 800,000 putative cells and 14.5 billion reads.

Workflow Overview

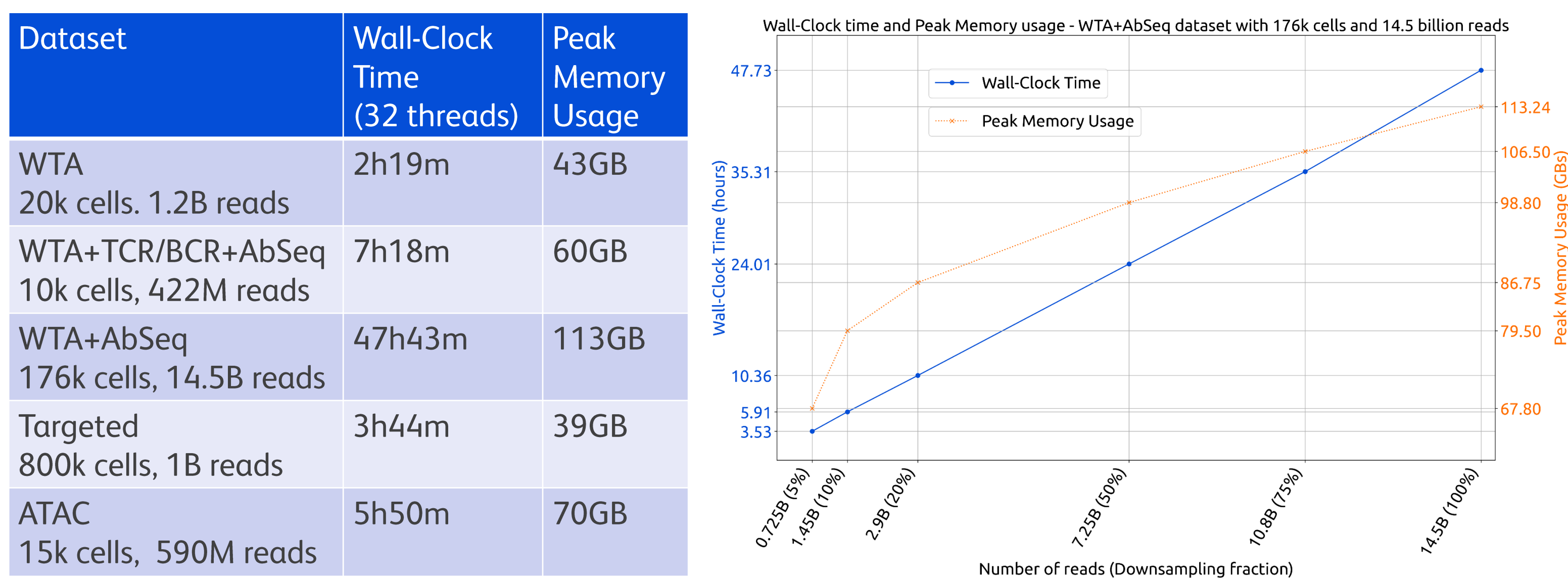


Performance Benchmarks

Comparison to previous pipeline release

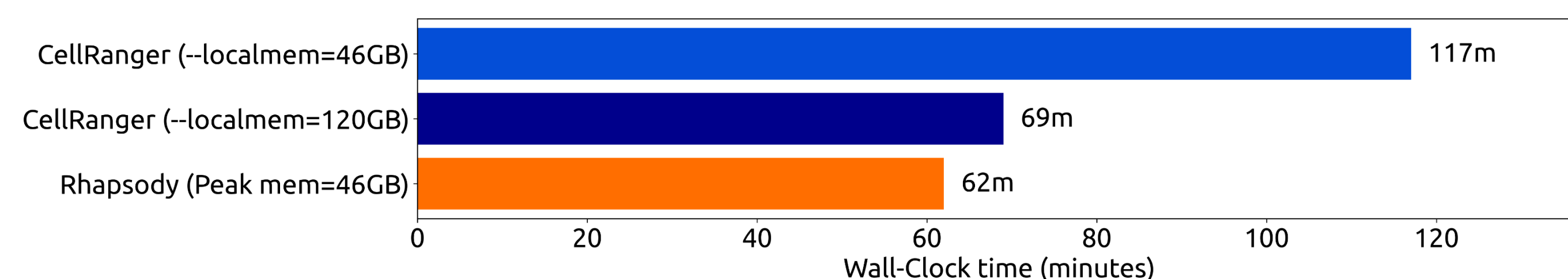


Runtime and Memory Usage



Note: All datasets were run on a m6i.8xlarge AWS instance with 32 CPUs, 128GB RAM and a 4 TB SSD EBS disk

Comparison b/w Rhapsody 2.2 and Cell Ranger 7.2.0 Dataset: 10k PBMCs, 496M WTA reads



Both programs were run with 32 threads

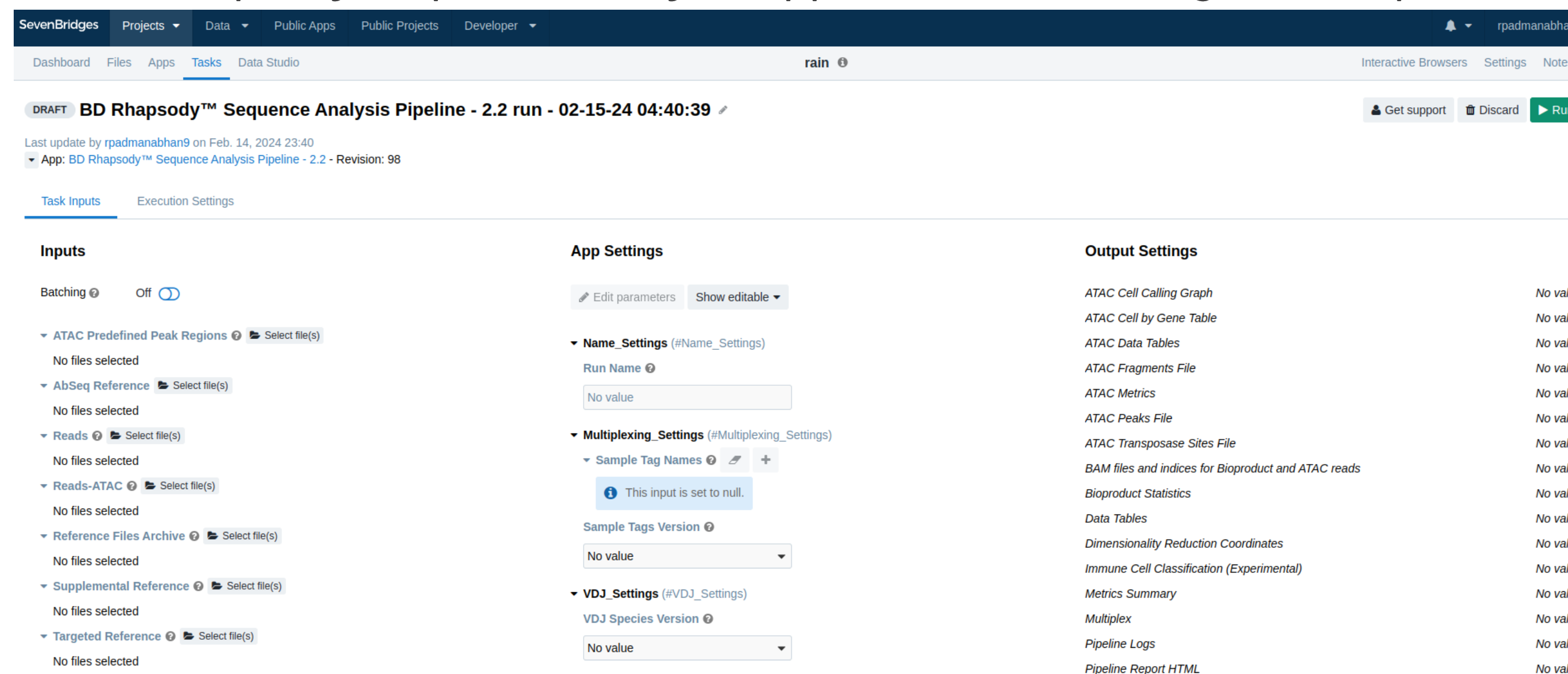
10X dataset: <https://www.10xgenomics.com/resources/datasets/10k-human-pbmc-3-v3-1-chromium-x-with-intronic-reads-3-1-high>

Rhapsody dataset: <https://bd-rhapsody-public.s3.amazonaws.com/Rhapsody-Demo-Data-Inputs/09WTA-ABC-EB-10kPBMC.tar> (09-WTA library downsampled to 496M reads)

Convenient deployment options

1. SevenBridges cloud platform
2. Local deployment using Docker container
3. Local deployment using Dockerfree bundle on Ubuntu 16.04+ / CentOS 8+ / RHEL 7+

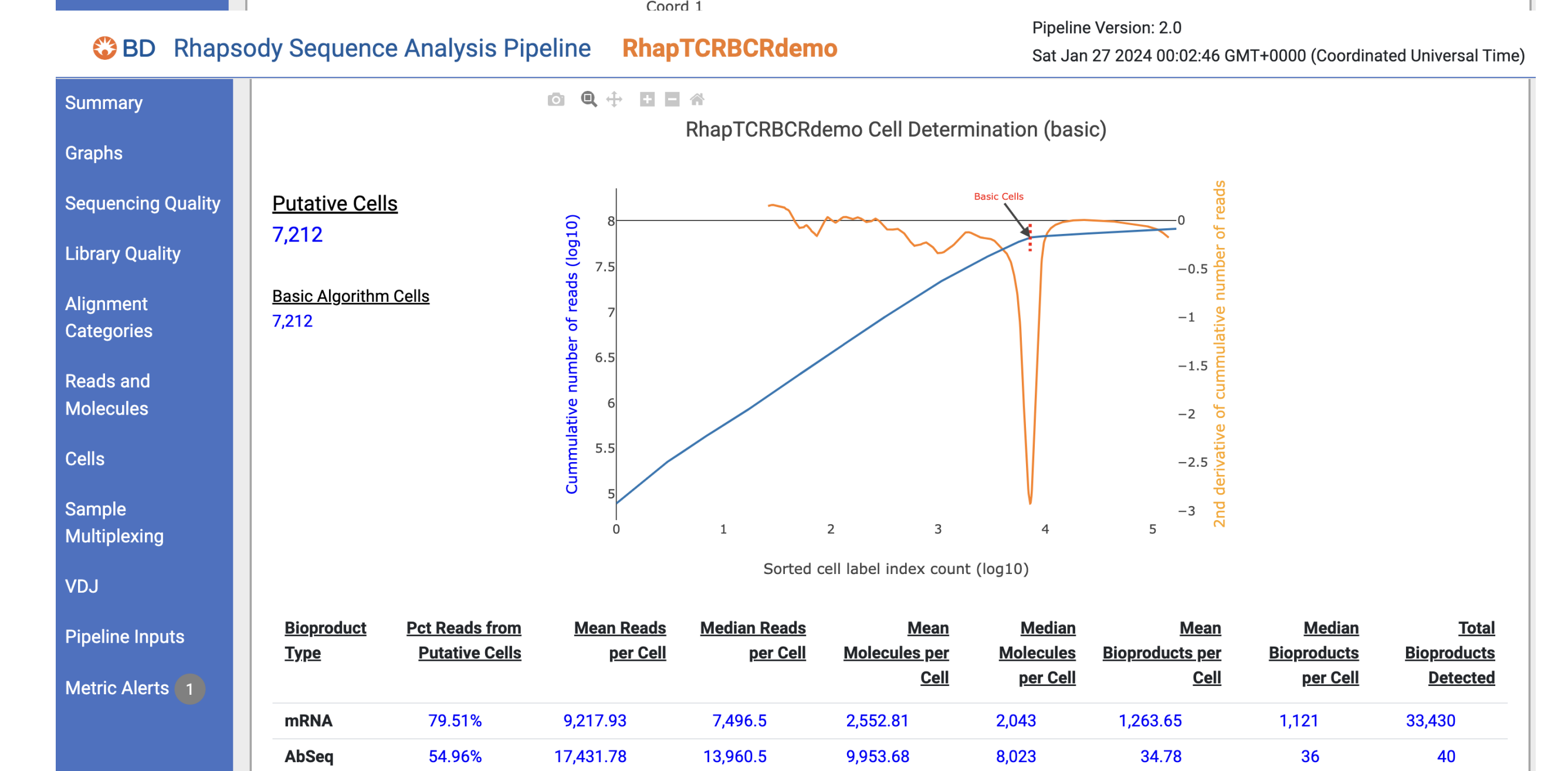
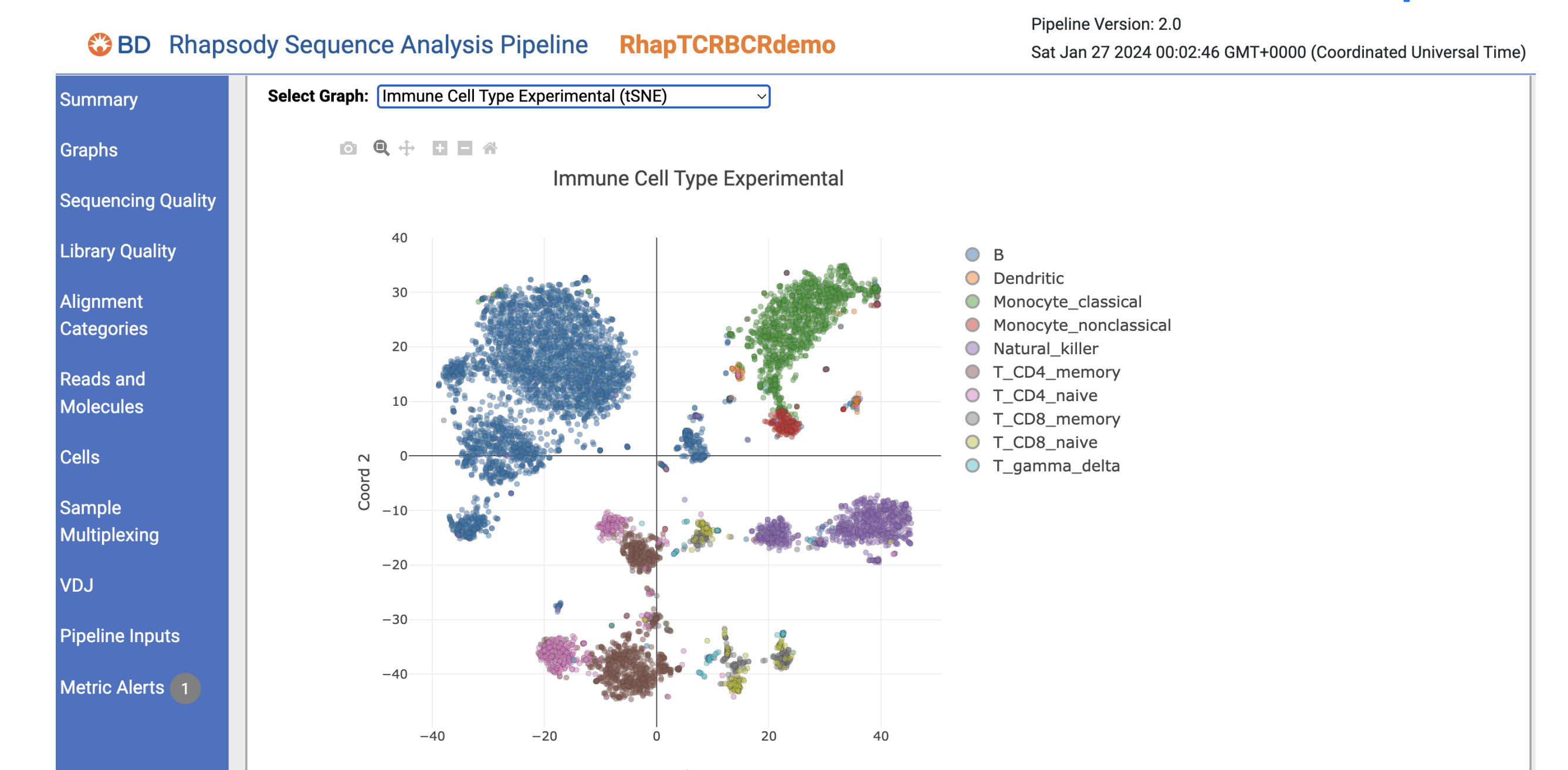
BD Rhapsody Sequence Analysis App on the SevenBridges cloud platform



Visit the link below to access the setup guide and documentation

<https://bd-rhapsody-bioinfo-docs.genomics.bd.com/>

Visualize results with the interactive HTML report



Output files are ready for downstream analysis

```
import muon
rhapsody_pbmc = muon.read_h5mu("Rhapsody-TCRBCR-Demo.h5mu")
rhapsody_pbmc

MuData object with n_obs x n_vars = 7212 x 33471
var: 'Raw_Reads', 'Raw_Molecules', 'Raw_Seq_Depth', 'RSEC_Adjusted_Molecules', 'RSEC_Adjusted_Reads_non-singleton', 'RSEC_2_modalities'
rna: 7212 x 33431
obs: 'Cell_Type_Experimental', 'Sample_Tag', 'Sample_Name', 'SampleTag01_hs_Read_Count', 'SampleTag02_hs_Read_Count'
var: 'Raw_Reads', 'Raw_Molecules', 'Raw_Seq_Depth', 'RSEC_Adjusted_Molecules', 'RSEC_Adjusted_Reads_non-singleton'
uns: 'Pipeline_Inputs', 'Pipeline_Metrics'
obsm: 'X_tsne', 'X_umap'
prot: 7212 x 40
obs: 'Cell_Type_Experimental', 'Sample_Tag', 'Sample_Name', 'SampleTag01_hs_Read_Count', 'SampleTag02_hs_Read_Count'
var: 'Raw_Reads', 'Raw_Molecules', 'Raw_Seq_Depth', 'RSEC_Adjusted_Molecules', 'RSEC_Adjusted_Reads_non-singleton'
uns: 'Pipeline_Inputs', 'Pipeline_Metrics'
obsm: 'X_tsne', 'X_umap'
```

For Research Use Only. Not for use in diagnostic or therapeutic procedures. BD, the BD Logo and BD Rhapsody are trademarks of Becton, Dickinson and Company or its affiliates. All other trademarks are the property of their respective owners. © 2024 BD. All rights reserved. BD-117497 (v1.0) 0324