

Introduction

The Genomics Research Group (GRG) generated data from hundreds of individual SUM149PT cells treated with the histone deacetylase inhibitor TSA vs. untreated controls across several scRNA-Seq platforms (Fluidigm C1, WaferGen iCell8, 10X Genomics Chromium Controller, and Illumina/BioRad ddSEQ). The goals of this project are to demonstrate RNA sequencing (RNA-Seq) methods for profiling the ultra-low amounts of RNA present in individual cells, and RNA amplification using the various currently available platforms. We will discuss the results of the study as well as technical challenges/lessons





sequencing reagents) and Illumina (Sequencing reagents, SureCell WTA reagents, ddSEQ demo, and dedicated support for SureCell analysis). Special thanks to Marcy Kuentzel at the Center for Functional Genomics at SUNY Albany and Michelle Zanche and Marlene Balys at the University of Rochester Medical Center for generating the samples. Thanks to Steven Lotz at the Neural Stem Cell Institute for assistance wit the iCell8 system. Thanks also to Baldwin Dilone at the Harvard Biopolymers facility for the single cell capture and library prep and Flow Cytometry Facility at URMC for cell sorting support.

Comparative analysis of single-cell RNA Sequencing Platforms and Methods

Sridar V. Chittur¹, Marcy Kuentzel¹, Michelle Zanche^{9,10}, Marlene Balys^{9,12}, Natalia Garcia-Reyero², Catharine Aquino Fournier³, Hubert Rehrauer³, Robert G. Steen⁴, Sarah A. Boswell⁴, Jin Jen⁵, Christopher E. Mason⁶, Jonathan Foox⁶, Shahin Mohammadi^{7,8}, Jose Davila-Velderrain^{7,8}, Manolis Kellis^{7,8}, Jason R. Myers^{9,10}, Jacquelyn Lillis^{9,10}, Andrew McDavid¹¹, John M Ashton^{9,10,12}

¹Center for Functional Genomics, University at Albany-SUNY; ²US Army ERDC Environmental Laboratory; ³Functional Genomics Center Zurich, ETHZ; ⁴Biopolymers Facility, Harvard Medical School; ⁵Genome Analysis Core, Mayo Clinic; ⁵Institute for Computational Biomedicine, Weill Cornell Medicine; ⁷Computer Science and Artifical Intelligence Laboratory (CSAIL), MIT, ⁸Broad Institute of MIT and Harvard; ⁹Department of Microbiology & Immunology, ¹⁰Genomics Research Center, University of Rochester, ¹¹Department of Biostatistics and Computational Biology, University of Rochester, ¹²James P. Wilmot **Cancer Institute, University of Rochester.**

Results & Observations

Platform	Technology	Capacity	Capture Efficiency	Cell Size Dependence	Cell Monitoring	UMI	Ease of Use	Type of Profiling	Cell Capture to Library Cost (per sample)	Cost (per Cell)	Cell Capture + Library Generation Time
erGen iCell8	Ordered Array	1,000-2,000	~30%	No	Yes	No	Moderate	3' expression	\$ 3,100.00	\$3.00	7hrs
digm 96	Nanofluidics	96	~20%	Yes	Yes	No	Moderate	whole transcript	\$ 2,000.00	\$20.83	10hrs
digm HT	Nanofluidics	400 - 800	~18%	Yes	Yes	No	Hard	3' expression	\$ 2,500.00	\$6.25 - 3.13	12hrs
Genomics v2	Droplet Encapsulation	1,000-10,000	~65%	No	No	Yes	Easy	3' expression	\$ 1,500.00	\$1.50 - \$0.15	10hrs
ina/BioRad ddSEQ	Droplet Encapsulation	300 - 1,200	~3%	No	No	Yes	Easy	3' expression	\$ 1,200.00	\$4.00 - \$1.00	10hrs

Table 1. Cost analysis and key attributes of scRNA-seq platforms. Costs were determined using list prices for each technology available through respective vendor.

Figure 2. Cell outlier detection using Scater automated parameters. Principal components for outlier detection results identify from Scater automated QC for data generated with Fluidigm C1 (A), WaferGen iCell8 (B), 10X Genomics with immediate processing (C), Fludigm HT (D), Illumina/BioRad ddSeq (E), or 10X Genomics with processing after overnight shipment (F). tSNE plots of single cell association based on treatment group (G).

Basic Metrics & Preliminary Assessment







Figure 3. TSA gene signature for Bulk RNA-Seq experimentation. PCA plot showing the relationship between DMSO and TSA treated SUM149PT cells from bulk TruSeq RNA-Seq (A). Volcano plot showing the extent of significant differentially expression genes between DMSO and TSA, p<0.05 (B). Heatmap of the 982 differentially expressed genes identified as statistically significant, p<0.05 (C).



Figure 4. Correlation among single-cell RNA-Seq platform technologies. Correlation of differential expression results between each platform and bulk RNA-Seq results (A). Correlation of differential expression results across platform technologies (B). Distribution of genes identified as differentially expression between DMSO vs TSA across all platforms tested (C).

 Each platform technology has some tradeoffs, such as in cell efficiency, data usability, or gene diversity.

Bulk RNA-Seq Differential Expression

Conclusions

• No platform reached gene detection saturation at sequencing depths tested, so all perform equally well in sensitivity under the tested conditions. • Some platforms showed higher degree of outliers and differing ability to resolve treatment groups well.

• Using bulk RNA-Seq (TruSeq) as "ground truth", a low concordance was observed for differentially expressed genes across platforms.