



Single-Cell ATAC-Seq

The SureCell® ATAC-Seq Library Prep Kit and ATAC-Seq Analysis Toolkit enable reproducible genome-wide profiling of the epigenomic landscape at the single-cell level with a high number of unique reads per cell so you can better understand the mechanisms that drive how genes are regulated.

The Bio-Rad single-cell ATAC-Seq (scATAC-Seq) solution provides users with:

- The ability to work with both whole cells and nuclei
- A high number of unique fragments that map to the nuclear genome, ATAC peaks, and transcription start sites (TSSs)
- Up to 90% cell capture efficiency
- A fast (<8 hr) simplified workflow with minimal hands-on time and convenient start and stop points
- Adjustable cell throughput range of as few as 400 or more than 4,000 cells per sample, more than 30,000 cells/kit
- A powerful and efficient informatics tool in a flexible pipeline

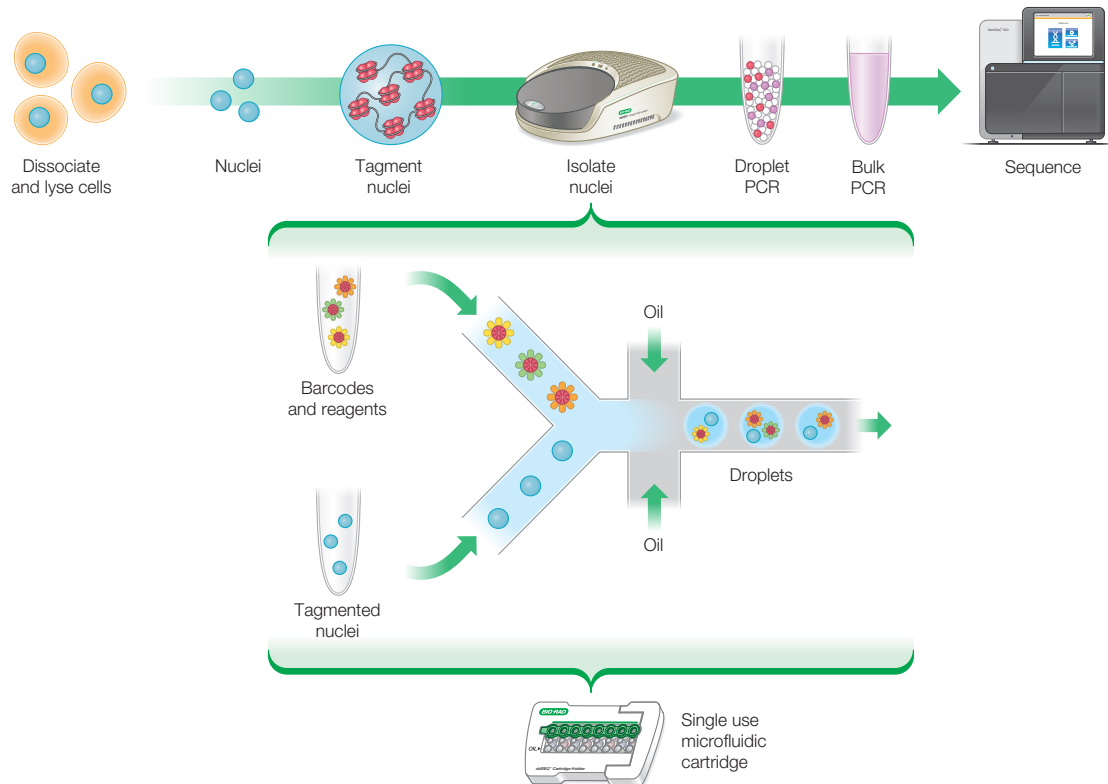


Visit [bio-rad.com/scATACSeqKit](https://www.bio-rad.com/scATACSeqKit) for more information.

High-Resolution Mapping of Open Chromatin with scATAC-Seq

Single-cell ATAC-Seq (assay for transposase-accessible chromatin using sequencing) with high cell throughput provides a novel method with which to map genome-wide chromatin accessibility for thousands of individual single cells. The simple one-day workflow provides easy access to epigenetic information, including transcriptional dynamics from cell to cell.

Single cells or their isolated nuclei are tagged at regions of open chromatin using hyperactive Tn5 transposase in a bulk reaction. Barcoding and amplification of tagged sites within individual droplets generate a library of fragments representative of the original open chromatin profile of each cell.



scATAC-Seq Workflow

Bio-Rad's scATAC-Seq workflow is streamlined and offers users:

- Less than 8 hours of cell and library prep time
- Sample-to-sample reproducibility
- Flexible control over workflow with multiple safe stop points
- Minimum hands-on time

1	Tagment isolated nuclei or whole cells	
2	Encapsulate tagmented nuclei/cells and barcode beads in droplets	
3	Barcode fragments	Safe Stop
4	Break emulsion and purify barcoded product	Safe Stop
5	Amplify barcoded fragments to generate libraries	

scATAC-Seq Applications

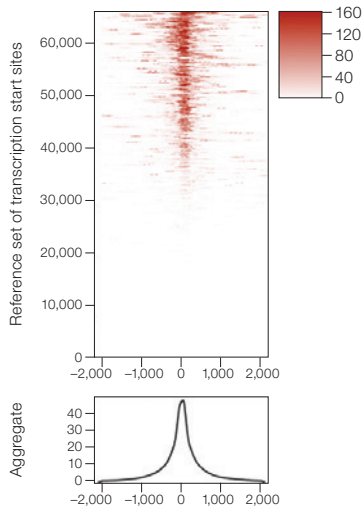
Single-cell ATAC-Seq allows you to take your studies to the next level.

Use scATAC-Seq to:

- Answer deep questions about regulation of gene expression through interrogation of promoters, enhancers, or other regulatory elements
- Validate single-cell RNA-Seq studies as an orthogonal approach
- Identify cell-specific regulatory signatures that are phenotypic of diseased cell variants and mechanisms of cell fate

Gene-Specific Data

High signal/noise with more than 40 times better alignment to TSSs



Bio-Rad's scATAC-Seq solution detects TSSs with high signal to noise. Reads are mapped to a reference set of TSSs to form an aggregate distribution aligned to TSSs and extended 2,000 bp in both directions. The value at the center of the distribution represents the TSS score.

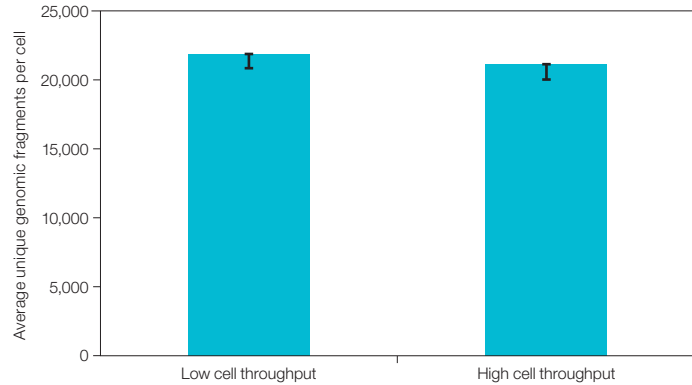
Capture Efficiency

Highly reproducible cell utilization can be obtained across multiple samples, experiments, or users.

Cell Input	Cell Output	Cell Utilization, %
7,392	4,644	63
7,840	5,073	65
7,488	4,621	62
7,648	4,443	58
6,848	4,340	63
6,656	4,649	70
6,976	5,044	72
6,400	3,720	58
6,976	4,351	62
7,456	3,735	50
3,184	3,159	99
6,380	4,477	70
6,022	3,777	63
7,424	5,116	69
6,976	4,331	62
7,424	4,786	64
7,193	4,240	59
8,128	6,130	75
8,384	6,506	78

Highly Sensitive and Reproducible Results

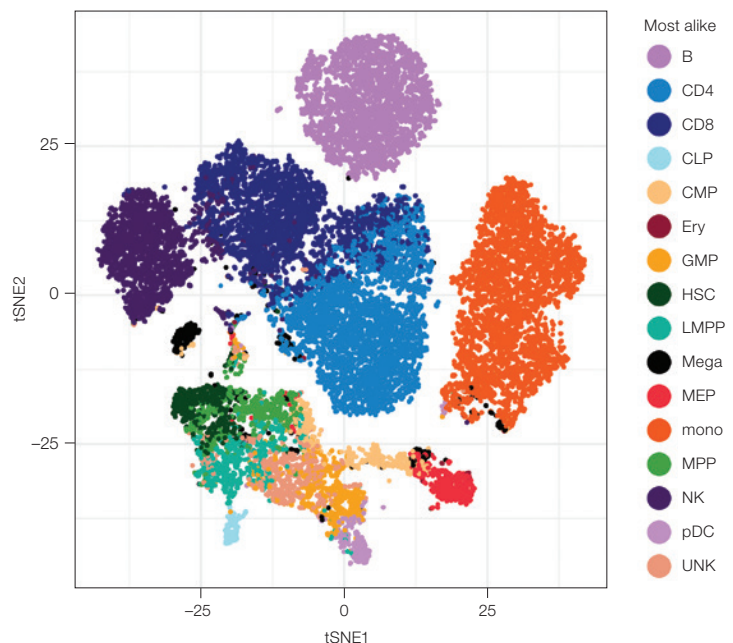
Achieve >20,000 unique fragments at 50,000 reads/cell for a wide dynamic range of cell throughput



	Mixed NIH 3T3 and K562 Cells	
	Low-Cell Throughput Samples	High-Cell Throughput Samples
Average unique genomic fragments per cell	21,828	21,176
Lower 95% confidence interval	20,761	19,956
	N = 42	N = 28

In a mixed-species experiment with an equal input of mouse NIH 3T3 and human K562 cells, greater than 20,000 unique genomic fragments per cell are detected at 50,000 reads/cell. Low-cell throughput samples had an average of 397 cells per sample, while high-cell throughput samples had an average of 4,408 cells per sample.

Cell Classification

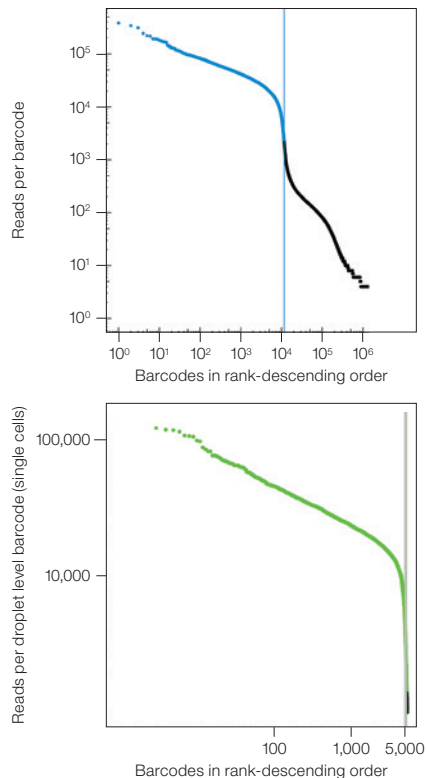


scATAC-Seq data can deconvolute highly complex cellular populations into their constitutive subsets. Using chromatin accessibility data, scATAC-Seq facilitates identifying and enumerating cell types within existing epigenomic data from complex cellular mixes. Analysis in collaboration with F. Duarte, C. Lareau, V. Kartha, Buenrostro Lab, Broad Institute of MIT and Harvard.

Simple, Yet Powerful, Data Analysis Options

Bio-Rad's bioinformatic pipeline enables rich data analytics in a transparent and flexible software pipeline of industry-standard tools. Secondary analysis starts with your FASTQ files and outputs biologically interpretable data that can be loaded into popular tertiary analysis tools like Seurat and ChromVAR.

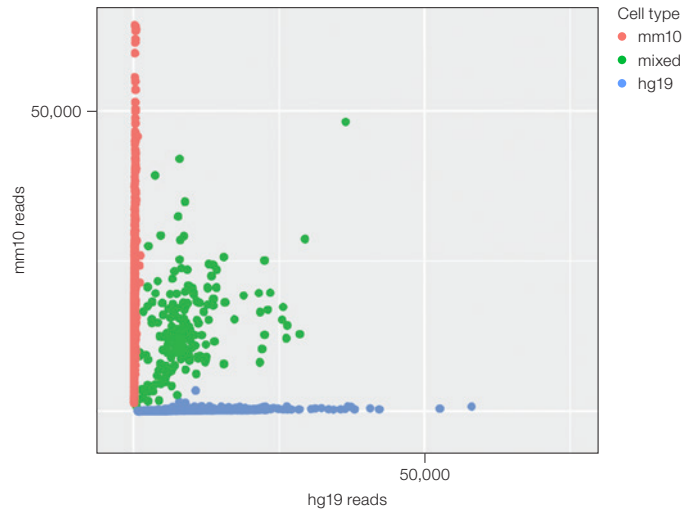
- Estimates gain and loss of chromatin accessibility within peaks
- Enables accurate and robust clustering of scATAC-Seq profiles and characterization of sequence motifs associated with gene expression
- Identifies cis- and trans-acting elements that are the source of diverse cellular phenotypes



Knee plots. Top, barcodes are sorted in descending order according to the number of unique reads per barcode. The data shown to the left of the vertical blue line represent barcodes that tag single cells. Pairwise comparisons are performed across these barcodes to identify those with a high percentage of shared fragments. These are merged bioinformatically to create droplet-level barcodes, which define single-cell data. Bottom, single cells sorted in descending order according to the number of unique reads per cell (green line).

Confident Single-Cell Transcript Identification

Minimum crosstalk (<5%) with more than 5,000 cells of mixed species



An equal number of mouse NIH 3T3 and human K562 cells were mixed and processed with the SureCell ATAC-Seq Library Prep Kit and ATAC-Seq Analysis Toolkit. These data show a sample with 5,024 cells. Droplets containing both human and mouse cells are shown in green and droplets with only human or mouse cells are shown in blue and red, respectively. Based on these data, cross talk, or the percentage of cell doublets, is calculated as 3.76%.

Ordering Information

Catalog #	Description
17004620	SureCell ATAC-Seq Library Prep Kit , includes SureCell ATAC-Seq Reagent Box A, SureCell ATAC-Seq Reagent Box B, and ddSEQ M Cartridges
12009359	ddSEQ M Cartridges
12009358	SureCell ATAC-Seq Reagent Box A
12009357	SureCell ATAC-Seq Reagent Box B
12009360	SureCell ATAC-Seq Index Kit
12004336	ddSEQ Single-Cell Isolator , includes instrument and associated component consumables

Visit bio-rad.com/scATACSeqKit for more information about single-cell ATAC-Seq and for additional application data.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc.

Purchase of Single-Cell NGS Sample Preparation products from Bio-Rad Laboratories includes a limited, non-transferable right to use the product only to perform internal research in the field of single-cell analysis and with Bio-Rad's ddSEQ Single-Cell Isolator for research purposes only. Unless we expressly state otherwise in additional Terms and Conditions, no rights are granted for you to distribute or resell the products or to use the products for commercial applications, such as contract services or fee for services. No rights are granted for diagnostic uses. Information concerning a license for such uses can be obtained from Bio-Rad Laboratories. It is the responsibility of the purchaser/end user to acquire any additional intellectual property rights that may be required. The Single-Cell NGS Sample Preparation products and ddSEQ System and/or their use is covered by claims of U.S. patents, and/or pending U.S. and non-U.S. patent applications owned by or under license to Bio-Rad Laboratories, Inc. See bio-rad.com/en-us/trademarks for details.

SureCell and the corresponding logos are trademarks of Illumina, Inc. in the U.S. and other countries, and are used under license. All trademarks used herein are the property of their respective owner.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550 **Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00 **Germany** 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30 **Poland** 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580 **Sweden** 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2576 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 01923 47 1301

