

Single-Cell ATAC-Seq

The ddSEQ™ 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:

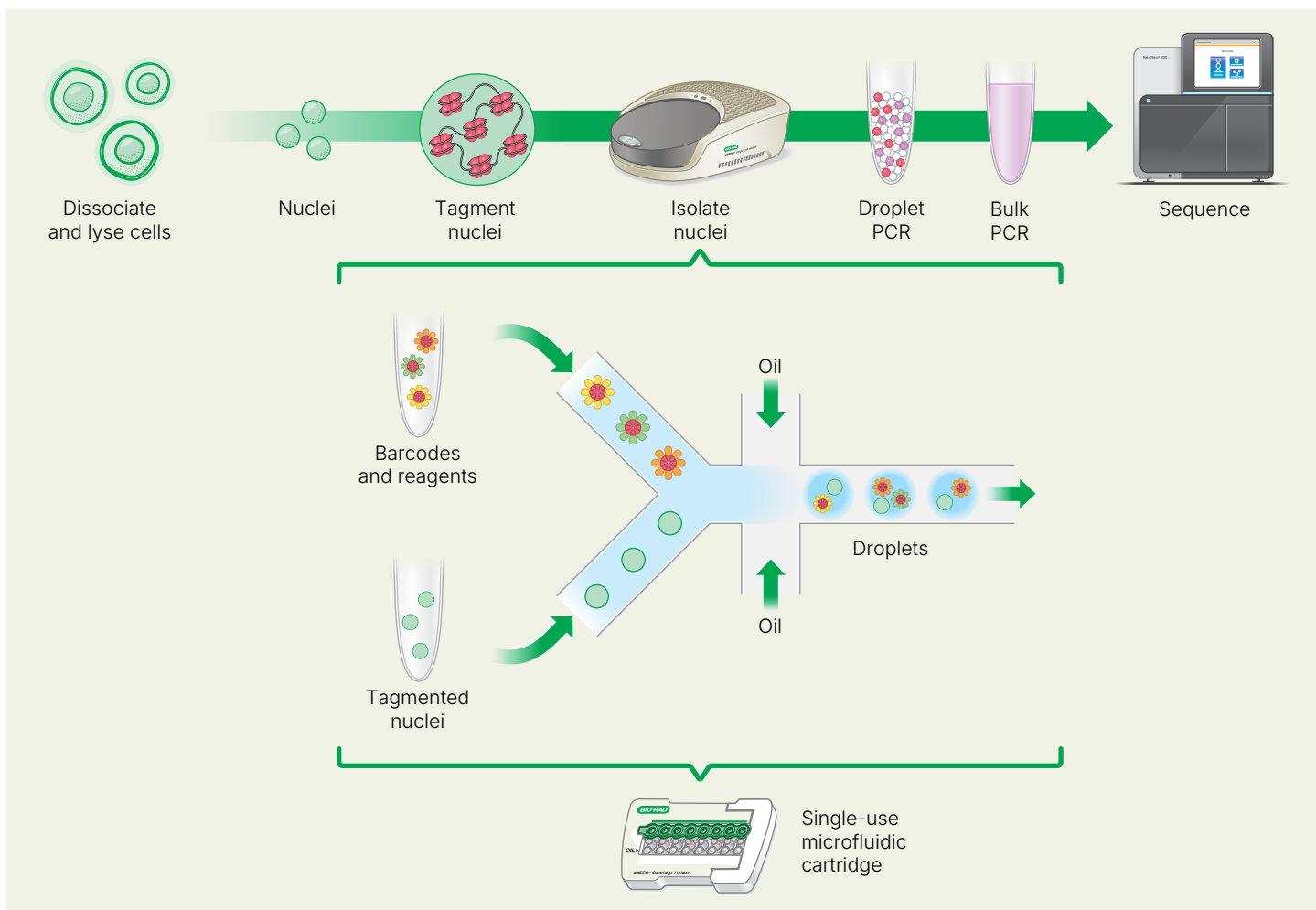
- The ability to work with both whole cells and nuclei
- A high number of unique fragments that map to the nuclear genome, ATAC peaks, and transcription start sites (TSSs)
- Up to 90% cell capture efficiency
- A fast (<8 hr) simplified workflow with minimal hands-on time and convenient start and stop 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



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

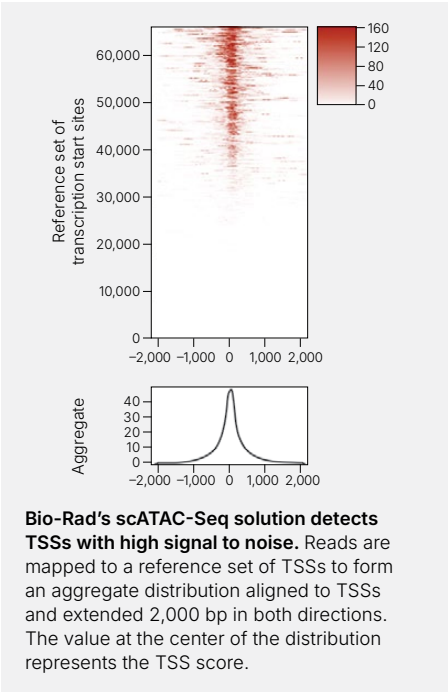
Single-cell ATAC-Seq allows you to take your studies to the next level.

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

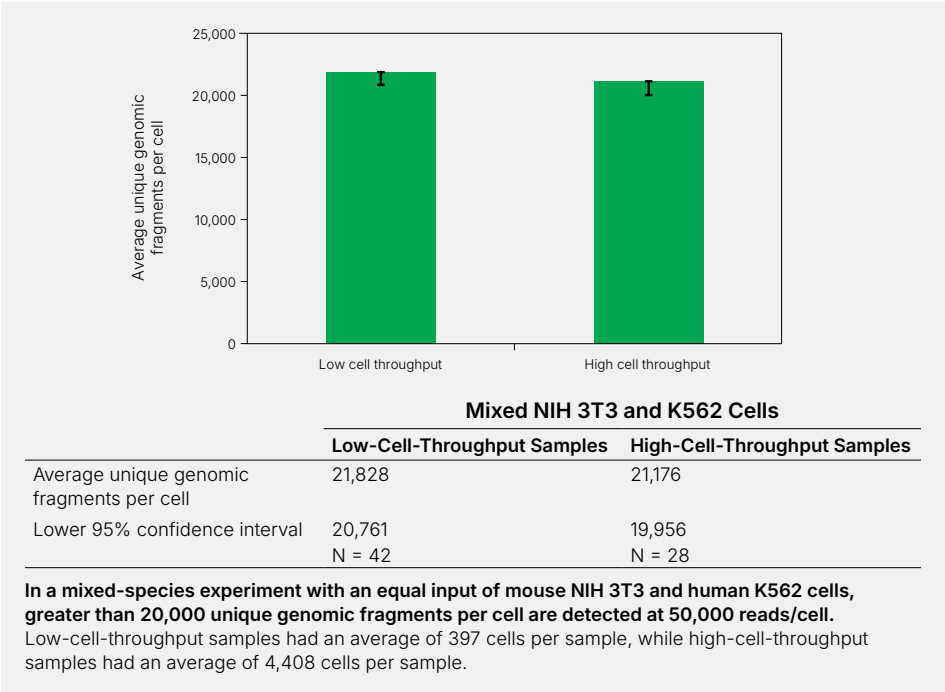
Gene-Specific Data

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



Highly Sensitive and Reproducible Results

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

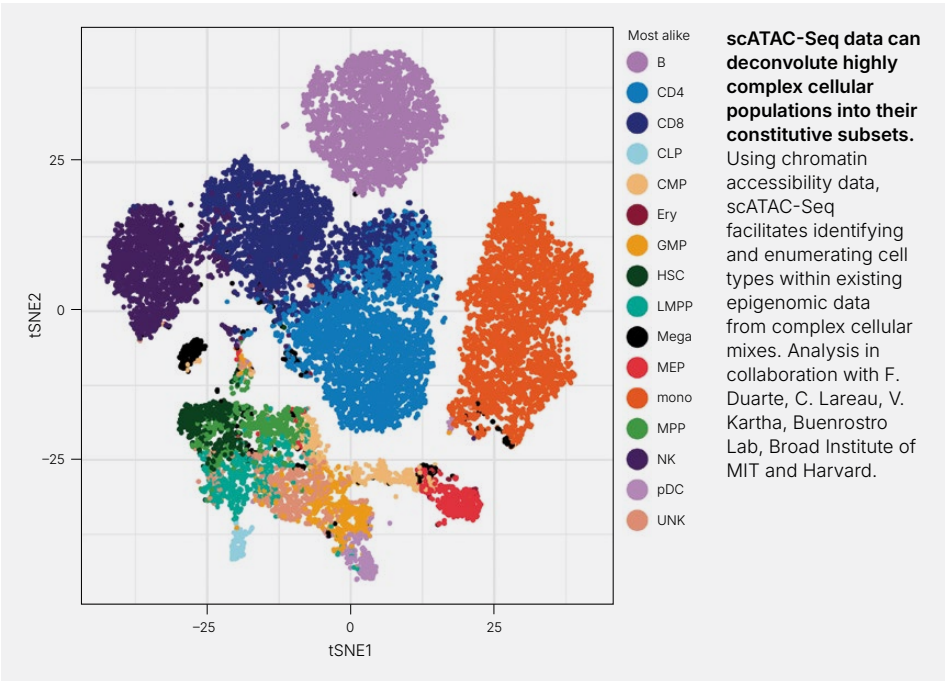


Capture Efficiency

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

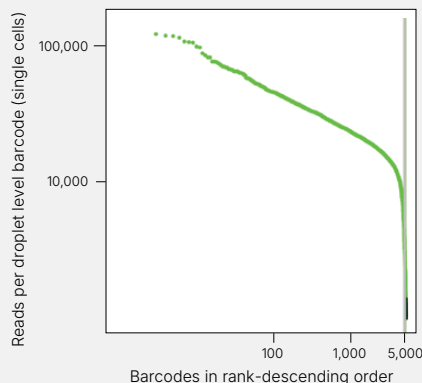
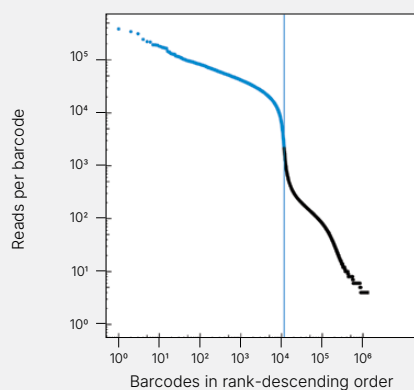
Cell Classification



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.

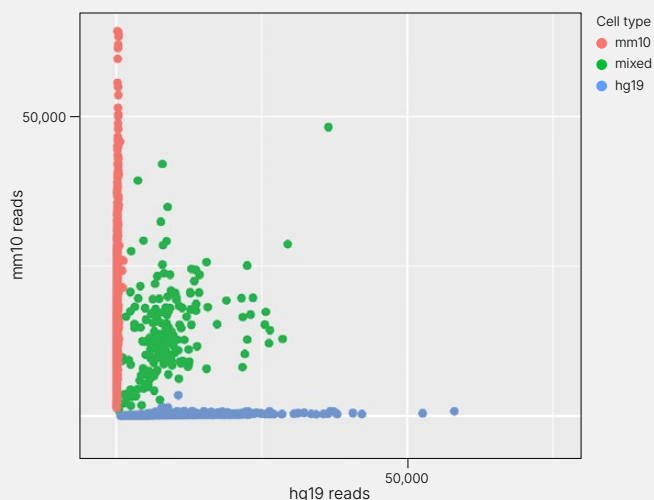
- Estimates gain and loss of chromatin accessibility within peaks
- Enables accurate and robust clustering of scATAC-Seq profiles and characterization of sequence motifs associated with gene expression
- Identifies cis- and trans-acting elements that are the source of diverse cellular phenotypes



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

Visit [bio-rad.com/scATACSeqKit](https://www.bio-rad.com/scATACSeqKit) for more information about single-cell ATAC-Seq and for additional application data.

Bio-Rad and ddSEQ are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.

© 2026 Bio-Rad Laboratories, Inc.

Purchase of Digital PCR and/or Single-Cell NGS Sample Preparation products (the "Products") from Bio-Rad Laboratories is subject to Bio-Rad Laboratories, Inc. Standard Terms and Conditions of Sale, which can be accessed at <https://www.bio-rad.com/en-us/terms-conditions>. Unless we expressly state otherwise in additional Terms and Conditions, no rights are granted for you to distribute or resell the Products. Unless we expressly state otherwise in additional Terms and Conditions, no rights are granted for the development or commercialization of diagnostic assays for use with the Products without a license from Bio-Rad. It is the user's obligation to obtain a commercial license from Bio-Rad for (i) all commercial uses (not just diagnostic uses) and (ii) sale of assays for use on Bio-Rad's dPCR and ddSEQ instruments. The Products and/or their use are covered by U.S. and foreign patents and/or pending patent applications owned by or under license to Bio-Rad Laboratories, Inc. See <https://www.bio-rad.com/en-us/trademarks>.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website [bio-rad.com](https://www.bio-rad.com) USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 4003 0399 Canada 1 800 361 1808 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23 France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 0 3 9636050 Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 080 007 7373 Luxembourg 00 800 00 24 67 23 Mexico 52 555 488 7670 The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Poland 00 800 00 24 67 23 Portugal 00 800 00 24 67 23 Russian Federation 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23 Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23 Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23