

Droplet Digital Sequencing (ddSEQ) Solution Publications List



ddSEQ Technology

Bulletin 7168

The ddSEQ workflow utilizes a droplet-based single-cell sequencing technology, enabling cancer, immunology, and stem cell research in various fields.

Different applications help researchers assess cell-to-cell heterogeneity, identify new subpopulations, and discover epigenetic markers. Below is a collection of research articles that highlight the advantages and applications of the ddSEQ Single-Cell Isolator.



Single-Cell RNA Sequencing

Bernard V et al. (2019).

Single-cell transcriptomics of pancreatic cancer precursors demonstrates epithelial and microenvironmental heterogeneity as an early event in neoplastic progression. Clin Cancer Res 25, 2,194–2,205.



Tran KA et al. (2019).

Defining reprogramming checkpoints from single-cell analyses of induced pluripotency. Cell Rep 27, 1,726–1,741.



Nguyen A et al. (2018).

Single cell RNA sequencing of rare immune cell populations. Front Immunol 9, 1,553.



Sarkar A et al. (2018).

Efficient generation of CA3 neurons from human pluripotent stem cells enables modeling of hippocampal connectivity in vitro. Cell Stem Cell 22, 684–697.



Shih AJ et al. (2018).

Identification of grade and origin specific cell populations in serous epithelial ovarian cancer by single cell RNA-seq. PLoS One 13, e0206785.





Single-Cell ATAC Sequencing

Lareau CA et al. (2019).

Droplet-based combinatorial indexing for massive-scale single-cell chromatin accessibility.

Nat Biotechnol 37, 916–924.



Visit [bio-rad.com/ddSEQ](https://www.bio-rad.com/ddSEQ) for more information about the ddSEQ solution.

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

All trademarks used herein are the property of their respective owner.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site [bio-rad.com](https://www.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 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 01923 47 1301

