



ZE5 Cell Analyzer

High Impact Publications List

Cancer Research

Au L et al. (2021).

Cytokine release syndrome in a patient with colorectal cancer after vaccination with BNT162b2.
Nat Med 27, 1,362–1,366.

Han J et al. (2021).

Resident and circulating memory T cells persist for years in melanoma patients with durable responses to immunotherapy.
Nat Cancer 2, 300–311.

Kieffer Y et al. (2020).

Single-cell analysis reveals fibroblast clusters linked to immunotherapy resistance in cancer.
Cancer Discov 10, 1,330–1,351.

Nam J et al. (2021).

Modularly programmable nanoparticle vaccine based on polyethyleneimine for personalized cancer immunotherapy.
Adv Sci (Weinh) 8, 2002577.

Perry JM et al. (2020).

Overcoming Wnt- β -catenin dependent anticancer therapy resistance in leukaemia stem cells.
Nat Cell Biol 22, 689–700.

Cell and Gene Therapy

Evgin L et al. (2022).

Oncolytic virus-mediated expansion of dual-specific CAR T cells improves efficacy against solid tumors in mice.
Sci Transl Med 14, eabn2231.

Kumar R et al. (2020).

Efficient polymer-mediated delivery of gene-editing ribonucleoprotein payloads through combinatorial design, parallelized experimentation, and machine learning.
ACS Nano [online ahead of print Nov 23, 2020]. Accessed August 17, 2022.

Lopez de Lapuente Portilla A et al. (2022).

Genome-wide association study on 13,167 individuals identifies regulators of blood CD34⁺ cell levels.
Blood 139, 1,659–1,669.

Mishra AK et al. (2021).

Preclinical development of CD126 CAR-T cells with broad antitumor activity.
Blood Cancer J 11, 3.

O'Neal J et al. (2022).

CS1 CAR-T targeting the distal domain of CS1 (SLAMF7) shows efficacy in high tumor burden myeloma model despite fratricide of CD8⁺CS1 expressing CAR-T cells.
Leukemia 36, 1,625–1,634.

Immunology

Ma S et al. (2021).

A role of PIEZO1 in iron metabolism in mice and humans.
Cell 184, 969–982.e13.

Nasrallah R et al. (2020).

A distal enhancer at risk locus 11q13.5 promotes suppression of colitis by T_{reg} cells.
Nature 583, 447–452.

Pasciuto E et al. (2020).

Microglia require CD4 T cells to complete the fetal-to-adult transition.
Cell 182, 625–640.e24.

Sawada Y et al. (2021).

Cutaneous innate immune tolerance is mediated by epigenetic control of MAP2K3 by HDAC8/9.
Sci Immunol 6, eabe1935.

Yeung F et al. (2020).

Altered immunity of laboratory mice in the natural environment is associated with fungal colonization.
Cell Host Microbe 27, 809–822.e6.

Infectious Disease

Chen RE et al. (2021).

Resistance of SARS-CoV-2 variants to neutralization by monoclonal and serum-derived polyclonal antibodies.
Nat Med 27, 717–726.

Grifoni A et al. (2020).

Targets of T cell responses to SARS-CoV-2 coronavirus in humans with COVID-19 disease and unexposed individuals.
Cell 181, 1,489–1,501.e15.

Lu Q et al. (2021).

SARS-CoV-2 exacerbates proinflammatory responses in myeloid cells through C-type lectin receptors and Tweety family member 2.
Immunity 54, 1,304–1,319.e9.

McCallum M et al. (2021).

N-terminal domain antigenic mapping reveals a site of vulnerability for SARS-CoV-2.
Cell 184, 2,332–2,347.e16.

Meng B et al. (2022).

Altered TMPRSS2 usage by SARS-CoV-2 Omicron impacts infectivity and fusogenicity.
Nature 603, 706–714.

Tarke A et al. (2022).

SARS-CoV-2 vaccination induces immunological T cell memory able to cross-recognize variants from Alpha to Omicron.
Cell 185, 847–859.e11.

Microbiology

Lin JD et al. (2020).

Rewilding Nod2 and Atg16l1 mutant mice uncovers genetic and environmental contributions to microbial responses and immune cell composition.
Cell Host Microbe 27, 830–840.e4.

Nakatsuji T et al. (2021).

Development of a human skin commensal microbe for bacteriotherapy of atopic dermatitis and use in a phase 1 randomized clinical trial.
Nat Med 27, 700–709.

Navaratna T et al. (2020).

Directed evolution using stabilized bacterial peptide display.
J Am Chem Soc 142, 1,882–1,894.

O'Neill AM et al. (2021).

Antimicrobials from a feline commensal bacterium inhibit skin infection by drug-resistant *S. pseudintermedius*.
Elife 10, e66793.

Tan R et al. (2022).

High-sugar, high-fat, and high-protein diets promote antibiotic resistance gene spreading in the mouse
intestinal microbiota.
Gut Microbes 14, 2022442.

Screening and Drug Discovery

Pinto D et al. (2020).

Cross-neutralization of SARS-CoV-2 by a human monoclonal SARS-CoV antibody.
Nature 583, 290–295.

Podracky CJ et al. (2021).

Laboratory evolution of a sortase enzyme that modifies amyloid- β protein.
Nat Chem Biol 17, 317–325.

Schardt JS et al. (2021).

Discovery and characterization of high-affinity, potent SARS-CoV-2 neutralizing antibodies via single B cell screening.
Sci Rep 11, 20738.

Starr TN et al. (2021).

SARS-CoV-2 RBD antibodies that maximize breadth and resistance to escape.
Nature 597, 97–102.

Tortorici MA et al. (2021).

Broad sarbecovirus neutralization by a human monoclonal antibody.
Nature 597, 103–108.

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

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner.
© 2022 Bio-Rad Laboratories, Inc.



**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 1 905 364 3435 **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 2 3473 4460 **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

