

Top 50 Journal Publications for the S3e Cell Sorter



Cell Sorting

Bulletin 7261



Cancer

Goossens S et al. (2017).

Oncogenic ZEB2 activation drives sensitivity toward KDM1A inhibition in T-cell acute lymphoblastic leukemia.
Blood 129, 981–990.



Logtenberg MEW et al. (2019).

Glutamyl cyclase is an enzymatic modifier of the CD47-SIRP α axis and a target for cancer immunotherapy.
Nat Med 25, 612–619.



Mezzadra R et al. (2017).

Identification of CMTM6 and CMTM4 as PD-L1 protein regulators.
Nature 549, 106–110.



Milanovic M et al. (2018).

Senescence-associated reprogramming promotes cancer stemness.
Nature 553, 96–100.



Montagner M et al. (2020).

Crosstalk with lung epithelial cells regulates Sfrp2-mediated latency in breast cancer dissemination.
Nat Cell Biol 22, 289–296.



Peirs S et al. (2017).

Targeting BET proteins improves the therapeutic efficacy of BCL-2 inhibition in T-cell acute lymphoblastic leukemia.
Leukemia 31, 2,037–2,047.



Cell Physiology

Rotty JD et al. (2017).

Arp2/3 complex is required for macrophage integrin functions but is dispensable for FcR phagocytosis and in vivo motility.
Dev Cell 42, 498–513.



Zhu H et al. (2017).

Impairments of spatial memory in an Alzheimer's disease model via degeneration of hippocampal cholinergic synapses.
Nat Commun 8, 1,676.





Genomics

Alexandrov A et al. (2017).

Fluorescence amplification method for forward genetic discovery of factors in human mRNA degradation.
Mol Cell 65, 191–201.

**Blomen VA et al. (2015).**

Gene essentiality and synthetic lethality in haploid human cells.
Science 350, 1,092–1,096.

**Brockmann M et al. (2017).**

Genetic wiring maps of single-cell protein states reveal an off-switch for GPCR signaling.
Nature 546, 307–311.

**Bulut-Karslioglu A et al. (2018).**

The transcriptionally permissive chromatin state of embryonic stem cells is acutely tuned to translational output.
Cell Stem Cell 22, 369–383.

**Gilsbach R et al. (2014).**

Dynamic DNA methylation orchestrates cardiomyocyte development, maturation and disease.
Nat Commun 5, 5,288.

**Loregger A et al. (2020).**

Haploid genetic screens identify SPRING/C12ORF49 as a determinant of SREBP signaling and cholesterol metabolism.
Nat Commun 11, 1,128.

**Ludwig AK et al. (2017).**

Binding of MBD proteins to DNA blocks Tet1 function thereby modulating transcriptional noise.
Nucleic Acids Res 45, 2,438–2,457.

**Manzo SG et al. (2018).**

DNA Topoisomerase I differentially modulates R-loops across the human genome.
Genome Biol 19, 100.

**Nothjunge S et al. (2017).**

DNA methylation signatures follow preformed chromatin compartments in cardiac myocytes.
Nat Commun 8, 1,667.

**Sheu-Gruttadauria J et al. (2019).**

Structural basis for target-directed microRNA degradation.
Mol Cell 75, 1,243–1,255.



**Tang WWC et al. (2015).**

A unique gene regulatory network resets the human germline epigenome for development.

Cell 161, 1,453–1,467.

**Vansant G et al. (2020).**

The chromatin landscape at the HIV-1 provirus integration site determines viral expression.

Nucleic Acids Res [published online ahead of print June 29, 2020].

Accessed July 27, 2020.



Immunology

Andrä I et al. (2020).

An evaluation of T-cell functionality after flow cytometry sorting revealed p38 MAPK activation.

Cytometry A 97, 171–183.

**Emgård J et al. (2018).**

Oxysterol sensing through the receptor GPR183 promotes the lymphoid-tissue-inducing function of innate lymphoid cells and colonic inflammation.

Immunity 48, 120–132.

**Gil-Cruz C et al. (2016).**

Fibroblastic reticular cells regulate intestinal inflammation via IL-15-mediated control of group 1 ILCs.

Nat Immunol 17, 1,388–1,396.

**Guo X et al. (2015).**

Innate lymphoid cells control early colonization resistance against intestinal pathogens through ID2-dependent regulation of the microbiota.

Immunity 42, 731–743.

**Lorentsen KJ et al. (2018).**

Bcl11b is essential for licensing Th2 differentiation during helminth infection and allergic asthma.

Nat Commun 9, 1,679.

**Morgantini C et al. (2019).**

Liver macrophages regulate systemic metabolism through non-inflammatory factors.

Nature Metabolism 1, 445–459.

**Pikor NB et al. (2020).**

Remodeling of light and dark zone follicular dendritic cells governs germinal center responses.

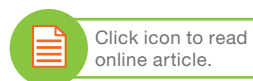
Nat Immunol 21, 649–659.

**Reisner K et al. (2017).**

Initiation of acute graft-versus-host disease by angiogenesis.

Blood 129, 2,021–2,032.





Microbiology

Duncan MC et al. (2019).

High-throughput analysis of gene function in the bacterial predator
Bdellovibrio bacteriovorus.
MBio 10, e01040-19.

**Duncan MC et al. (2018).**

Vibrio cholerae motility exerts drag force to impede attack by the bacterial predator
Bdellovibrio bacteriovorus.
Nat Commun 9, 4,757.

**Guiziu S et al. (2016).**

A part toolbox to tune genetic expression in *Bacillus subtilis*.
Nucleic Acids Res 44, 7,495–7,508.

**Huang L et al. (2018).**

Growth of *Mycobacterium tuberculosis* in vivo segregates with host macrophage
metabolism and ontogeny.
J Exp Med 215, 1,135–1,152.

**Kanjee U et al. (2017).**

CRISPR/Cas9 knockouts reveal genetic interaction between strain-transcendent
erythrocyte determinants of *Plasmodium falciparum* invasion.
Proc Natl Acad Sci USA 114, E9,356–E9,365.

**Lander N et al. (2015).**

CRISPR/Cas9-induced disruption of paraflagellar rod protein 1 and 2 genes in
Trypanosoma cruzi reveals their role in flagellar attachment.
MBio 6, e01012.

**Vilchèze C et al. (2017).**

Enhanced respiration prevents drug tolerance and drug resistance in
Mycobacterium tuberculosis.
Proc Natl Acad Sci USA 114, 4,495–4,500.

**Wu K et al. (2018).**

Role of PDGF receptor- α during human cytomegalovirus entry into fibroblasts.
Proc Natl Acad Sci USA 115, E9,889–E9,898.





Protein Analysis

Hill BJ et al. (2017).

A coiled-coil strategy for the directional display of multiple proteins on the surface of iron oxide nanoparticles.
RSC Adv 7, 12,133–12,143.

**Martínez-Limón A et al. (2016).**

Recognition of enzymes lacking bound cofactor by protein quality control.
Proc Natl Acad Sci USA 113, 12,156–12,161.

**Staus DP et al. (2020).**

Structure of the M2 muscarinic receptor- β -arrestin complex in a lipid nanodisc.
Nature 579, 297–302.

**Taylor ND et al. (2016).**

Engineering an allosteric transcription factor to respond to new ligands.
Nat Methods 13, 177–183.



Stem Cells

Borot F et al. (2019).

Gene-edited stem cells enable CD33-directed immune therapy for myeloid malignancies.
Proc Natl Acad Sci USA 116, 11,978–11,987.

**Chal J et al. (2015).**

Differentiation of pluripotent stem cells to muscle fiber to model Duchenne muscular dystrophy.
Nat Biotechnol 33, 962–969.

**Chen YW et al. (2017).**

A three-dimensional model of human lung development and disease from pluripotent stem cells.
Nat Cell Biol 19, 542–549.

**Cheng HW et al. (2019).**

Origin and differentiation trajectories of fibroblastic reticular cells in the splenic white pulp.
Nat Commun 10, 1,739.

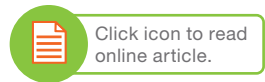
**Diaz-Cuadros M et al. (2020).**

In vitro characterization of the human segmentation clock.
Nature 580, 113–118.

**Irie N et al. (2015).**

SOX17 is a critical specifier of human primordial germ cell fate.
Cell 160, 253–268.





Kuroda T et al. (2019).

SALL3 expression balance underlies lineage biases in human induced pluripotent stem cell differentiation.
 Nat Commun 15, 2,175.



Murakami K et al. (2016).

NANOG alone induces germ cells in primed epiblast in vitro by activation of enhancers.
 Nature 529, 403–407.



Oberst P et al. (2019).

Temporal plasticity of apical progenitors in the developing mouse neocortex.
 Nature 573, 370–374.



Stuart HT et al. (2019).

Distinct molecular trajectories converge to induce naive pluripotency.
 Cell Stem Cell 25, 388–406.



Visit bio-rad.com/S3e for a wide range of videos, technical reports, published findings, and other resources describing Bio-Rad's S3e Cell Sorter.

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

Website 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

