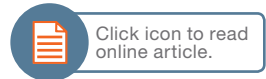


# Flow Cytometry Publications List



S3e™ and S3™ Cell Sorters

Bulletin 6741



Click icon to read online article.

## Cancer

Catenacci DV et al. (2015).

**Acquisition of portal venous circulating tumor cells from patients with pancreaticobiliary cancers by endoscopic ultrasound.**

Gastroenterology 149, 1,794–1,803.



Frame FM et al. (2016).

**Harvesting human prostate tissue material and culturing primary prostate epithelial cells.**

Methods Mol Biol 1443, 181–201.



Hernandez JR et al. (2015).

**Alternative CD44 splicing identifies epithelial prostate cancer cells from the mesenchymal counterparts.**

Med Oncol 32, 159.



Jin HJ et al. (2016).

**Identification and validation of regulatory SNPs that modulate transcription factor chromatin binding and gene expression in prostate cancer.**

Oncotarget 7, 54,616–54,626.



Kawahara T et al. (2016).

**Silodosin inhibits prostate cancer cell growth via ELK1 inactivation and enhances the cytotoxic activity of gemcitabine.**

Prostate 76, 744–756.



Lacotte S et al. (2016).

**Impact of myeloid-derived suppressor cell on Kupffer cells from mouse livers with hepatocellular carcinoma.**

Oncoimmunology 5, e1234565.



Li Y et al. (2015).

**P16INK4A is required for cisplatin resistance in cervical carcinoma SiHa cells.**

Oncol Lett 9, 1,104–1,108.





## Cancer *continued*

Peirs S et al. (2014).

**ABT-199 mediated inhibition of BCL-2 as a novel therapeutic strategy in T-cell acute lymphoblastic leukemia.**

Blood 124, 3,738–3,747.



Rae DT et al. (2015).

**A novel retroviral mutagenesis screen identifies prognostic genes in RUNX1 mediated myeloid leukemogenesis.**

Oncotarget 6, 30,664–30,674.



Rihani A et al. (2015).

**Inhibition of CDK4/6 as a novel therapeutic option for neuroblastoma.**

Cancer Cell Int 15, 76.



## Cell Characterization

Bezzerides VJ et al. (2016).

**CITED4 induces physiologic hypertrophy and promotes functional recovery after ischemic injury.**

JCI Insight 1, e85904.



Blomen VA et al. (2015).

**Gene essentiality and synthetic lethality in haploid human cells.**

Science 350, 1,092–1,096.



Hervier B et al. (2016).

**Involvement of NK cells and NKp30 pathway in antisynthetase syndrome.**

J Immunol 197, 1,621–1,630.



Konda N et al. (2016).

**Truncated EphA2 likely potentiates cell adhesion via integrins as well as infiltration and/or lodgment of a monocyte/macrophage cell line in the red pulp and marginal zone of the mouse spleen, where ephrin-A1 is prominently expressed in the vasculature.**

Histochem Cell Biol [published ahead of print September 24, 2016]. Accessed February 14, 2017.



Lothar A et al. (2016).

**Deoxycorticosterone acetate/salt-induced cardiac but not renal injury is mediated by endothelial mineralocorticoid receptors independently from blood pressure.**

Hypertension 67, 130–138.



Nührenberg TG et al. (2015).

**Cardiac myocyte de novo DNA methyltransferases 3a/3b are dispensable for cardiac function and remodeling after chronic pressure overload in mice.**

PLoS One 10, e0131019.





## Cell Characterization *continued*

Regn M et al. (2016).

**Peptidase inhibitor 16 is a membrane-tethered regulator of chemerin processing in the myocardium.**

J Mol Cell Cardiol 99, 57–64.



Turqueti-Neves A et al. (2015).

**The extracellular domains of IgG1 and T cell-derived IL-4/IL-13 are critical for the polyclonal memory IgE response in vivo.**

PLoS Biol 13, e1002290.



## Cell Development

Džinić T et al. (2016).

**Oxygen and differentiation status modulate the effect of X-ray irradiation on physiology and mitochondrial proteome of human neuroblastoma cells.**

Arch Physiol Biochem 122, 257–265.



Gilsbach R et al. (2014).

**Dynamic DNA methylation orchestrates cardiomyocyte development, maturation and disease.**

Nat Commun 5, 5,288.



Püngel S et al. (2015).

**Reconciling pillars of transient gene expression: From DNA prep via media, reagent and cell line development to holistic process optimization.**

BMC Proc 9, P18.



Saeki N et al. (2015).

**EphA2 promotes cell adhesion and spreading of monocyte and monocyte/macrophage cell lines on integrin ligand-coated surfaces.**

Cell Adh Migr 9, 469–482.



## Cell Differentiation

Afzal E et al. (2013).

**Nanolipodendrosome-loaded glatiramer acetate and myogenic differentiation 1 as augmentation therapeutic strategy approaches in muscular dystrophy.**

Int J Nanomedicine 8, 2,943–2,960.



Freedman BD et al. (2013).

**Adrenocortical zonation results from lineage conversion of differentiated zona glomerulosa cells.**

Dev Cell 26, 666–673.





## Cell Differentiation *continued*

Nam S et al. (2016).

**Interferon regulatory factor 4 (IRF4) controls myeloid-derived suppressor cell (MDSC) differentiation and function.**

J Leukoc Biol 100, 1,273–1,284.



Onder L et al. (2015).

**Alternative NF- $\kappa$ B signaling regulates mTEC differentiation from podoplanin-expressing precursors in the cortico-medullary junction.**

Eur J Immunol 45, 2,218–2,231.



Sécca C et al. (2016).

**IRF2BP2 transcriptional repressor restrains naive CD4 T cell activation and clonal expansion induced by TCR triggering.**

J Leukoc Biol 100, 1,081–1,091.



## Cell Identification

Carulli G et al. (2015).

**Combination of CD157 and FLAER to detect peripheral blood eosinophils by multiparameter flow cytometry.**

J Clin Exp Hematop 55, 55–60.



## Cell Sorting

Riddell A et al. (2015).

**Rmax: A systematic approach to evaluate instrument sort performance using center stream catch.**

Methods 82, 64–73.



## Epigenetics

Preissl S et al. (2015).

**Deciphering the epigenetic code of cardiac myocyte transcription.**

Circ Res 117, 413–423.



Tang WW et al. (2015).

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

Cell 161, 1,453–1,467.





## Immunology

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.



Imhof BA et al. (2016).

**CCN1/CYR61-mediated meticulous patrolling by Ly6Clow monocytes fuels vascular inflammation.**

Proc Natl Acad Sci USA 113, E4,847–E4,856.



Menendez CM et al. (2016).

**Resident T cells are unable to control herpes simplex virus-1 activity in the brain ependymal region during latency.**

J Immunol 197, 1,262–1,275.



## Microbial Analysis

Bidlingmaier S et al. (2016).

**Proteome-wide identification of novel ceramide-binding proteins by yeast surface cDNA display and deep sequencing.**

Mol Cell Proteomics 15, 1,232–1,245.



Eshaghi M et al. (2016).

**Brighter fluorescent derivatives of UTI89 utilizing a monomeric vGFP.**

Pathogens 5, E3.



Guiziou S et al. (2016).

**A part toolbox to tune genetic expression in *Bacillus subtilis*.**

Nucleic Acids Res 44, 7,495–7,508.



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.



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–e01015.



Liu Y et al. (2016).

**Immune activation of the host cell induces drug tolerance in *Mycobacterium tuberculosis* both in vitro and in vivo.**

J Exp Med 213, 809–825.



Moore SJ et al. (2016).

**EcoFlex: A multifunctional MoClo kit for *E. coli* synthetic biology.**

ACS Synth Biol 5, 1,059–1,069.





## Protein Analysis

Eshaghi M et al. (2015).

**Rational structure-based design of bright GFP-based complexes with tunable dimerization.**

Angew Chem Int Ed Engl 54, 13,952–13,956.



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.



Whiten DR et al. (2016).

**Rapid flow cytometric measurement of protein inclusions and nuclear trafficking.**

Sci Rep 6, 31138.



## Stem Cells

Bertacchi M et al. (2015).

**Activin/nodal signaling supports retinal progenitor specification in a narrow time window during pluripotent stem cell neuralization.**

Stem Cell Reports 5, 532–545.



Chal J et al. (2015).

**Differentiation of pluripotent stem cells to muscle fiber to model Duchenne muscular dystrophy.**

Nat Biotechnol 33, 962–969.



Irie N and Surani MA (2017).

**Efficient induction and isolation of human primordial germ cell-like cells from competent human pluripotent stem cells.**

Methods Mol Biol 1463, 217–226.



Murakami K et al. (2016).

**NANOG alone induces germ cells in primed epiblast in vitro by activation of enhancers.**

Nature 529, 403–407.



Pacini S et al. (2016).

**Mesangiogenic progenitor cells derived from one novel CD64(bright) CD31(bright)CD14(neg) population in human adult bone marrow.**

Stem Cells Dev 25, 661–673.



Quintanilla RH Jr et al. (2016).

**Kinetic measurement and real time visualization of somatic reprogramming.**

J Vis Exp, issue 113 [video].





## Virology

Browning DL et al. (2016).

**Insulated foamy viral vectors.**

Hum Gene Ther 27, 255–266.



Ramanujam D et al. (2016).

**Viral vector-based targeting of miR-21 in cardiac nonmyocyte cells reduces pathologic remodeling of the heart.**

Mol Ther 24, 1,939–1,948.



Yadav J et al. (2014).

**Opportunistic infections and complications in human immunodeficiency virus-1-infected children: Correlation with immune status.**

Sultan Qaboos Univ Med J 14, e513–e521.



## Flow Cytometry Posters

Lakshmiopathy U et al.

**Real time visualization and kinetic measurement of somatic reprogramming.**

Presented at: International Society for Stem Cell Research; June 24–27, 2015; Stockholm, Sweden.



Litterst L et al.

**Rapid and ultra-sensitive single-cell transcript profiling with Droplet Digital PCR (ddPCR™): Application to neuronal differentiation.**

Presented at: Neuroscience 2014; November 15–19, 2014; Washington, D.C.



Riddell A.

**Using a spectra physics Tsunami infrared pulsed laser and second harmonic generation on a Beckman Coulter MoFlo Legacy sorter: A tunable light source.**

Presented at: 30th Congress of the International Society for Advancement of Cytometry; June 26–30, 2015; Glasgow, Scotland. Abstract 421/B290, 248.



Riddell A et al.

**Sorting and maintaining haploid stem cells using a Beckman Coulter MoFlo Legacy and a Bio-Rad S3 by light scatter.**

Presented at: 30th Congress of the International Society for Advancement of Cytometry; June 26–30, 2015; Glasgow, Scotland. Abstract 420/B289, 247–248.



Visit [bio-rad.com/web/S3ePublished](http://bio-rad.com/web/S3ePublished) for a wide range of videos, technical reports, published findings, and other resources describing Bio-Rad's S3e Cell Sorter.



**Bio-Rad  
Laboratories, Inc.**

*Life Science  
Group*

**Web site** [bio-rad.com](http://bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 177 **Belgium** 32 (0)3 710 53 00 **Brazil** 55 11 3065 7550  
**Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 45 44 52 10 00 **Finland** 358 09 804 22 00  
**France** 33 01 47 95 69 65 **Germany** 49 89 31 884 0 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300  
**Israel** 972 03 963 6050 **Italy** 39 02 216091 **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 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04  
**Singapore** 65 6415 3188 **South Africa** 27 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 46 08 555 12700 **Switzerland** 41 026674 55 05  
**Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 020 8328 2000

