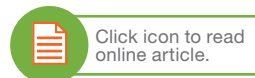


# Publication List: Electroporation of Nucleic Acids and Proteins for CRISPR-Related Gene Editing



Gene Pulser® / Gene Pulser Xcell™ / MicroPulser™ Electroporation Systems

Bulletin 6826



Click icon to read online article.

## 2016

### Highly efficient mouse genome editing by CRISPR ribonucleoprotein electroporation of zygotes.

Chen S, Lee B, Lee AY, Modzelewski AJ, He L (2016). J Biol Chem [published online ahead of print May 5, 2016]. Accessed June 14, 2016



### Chemical control of grafted human PSC-derived neurons in a mouse model of Parkinson's disease.

Chen Y, Xiong M, Dong Y, Haberman A, Cao J, Liu H, Zhou W, Zhang SC (2016). Cell Stem Cell 18, 817–826.



### Combinatorial metabolic pathway assembly in the yeast genome with RNA-guided Cas9.

EauClaire SF, Zhang J, Rivera CG, Huang LL (2016). J Ind Microbiol Biotechnol 43, 1,001–1,015.



### Human monocytes engage an alternative inflammasome pathway.

Gaidt MM, Ebert TS, Chauhan D, Schmidt T, Schmid-Burgk JL, Rapino F, Robertson AA, Cooper MA, Graf T, Hornung V (2016). Immunity 44, 833–846.



### Activation of RNase L is dependent on OAS3 expression during infection with diverse human viruses.

Li Y, Banerjee S, Wang Y, Goldstein SA, Dong B, Gaughan C, Silverman RH, Weiss SR (2016). Proc Natl Acad Sci USA 113, 2,241–2,246.



### Smad2 and Smad3 have differential sensitivity in relaying TGFβ signaling and inversely regulate early lineage specification.

Liu L, Liu X, Ren X, Tian Y, Chen Z, Xu X, Du Y, Jiang C, Fang Y, Liu Z, Fan B, Zhang Q, Jin G, Yang X, Zhang X (2016). Sci Rep 6, 21602.



### The CRISPR RNA-guided surveillance complex in *Escherichia coli* accommodates extended RNA spacers.

Luo ML, Jackson RN, Denny SR, Tokmina-Lukaszewska M, Maksimchuk KR, Lin W, Bothner B, Wiedenheft B, Beisel CL (2016). Nucleic Acids Res [published online ahead of print May 12, 2016]. Accessed June 14, 2016.



### Pathogenic variants in PIGG cause intellectual disability with seizures and hypotonia.

Makrythanasis P, Kato M, Zaki MS, Saito H, Nakamura K, Santoni FA, Miyatake S, Nakashima M, Issa MY, Guipponi M, Letourneau A, Logan CV, Roberts N, Parry DA, Johnson CA, Matsumoto N, Hamamy H, Sheridan E, Kinoshita T, Antonarakis SE, Murakami Y (2016). Am J Hum Genet 98, 615–626.



---

**Alternative splicing of MALT1 controls signalling and activation of CD4+ T cells.**

Meininger I, Griesbach RA, Hu D, Gehring T, Seeholzer T, Bertossi A, Kranich J, Oeckinghaus A, Eitelhuber AC, Greczmiel U, Gewies A, Schmidt-Supprian M, Ruland J, Brocker T, Heissmeyer V, Heyd F, Krappmann D (2016). Nat Commun 7, 11292.



---

**The SAGA deubiquitination module promotes DNA repair and class switch recombination through ATM and DNAPK-mediated  $\gamma$ H2AX formation.**

Ramachandran S, Haddad D, Li C, Le MX, Ling AK, So CC, Nepal RM, Gommerman JL, Yu K, Ketela T, Moffat J, Martin A (2016). Cell Rep 15, 1,554–1,565.



---

**A non-inheritable maternal Cas9-based multiple-gene editing system in mice.**

Sakurai T, Kamiyoshi A, Kawate H, Mori C, Watanabe S, Tanaka M, Uetake R, Sato M, Shindo T (2016). Sci Rep 6, 20011.



---

**Knockdown of *EPHA1* by CRISPR/CAS9 promotes adhesion and motility of HRT18 colorectal carcinoma cells.**

Wu BO, Jiang WG, Zhou D, Cui YX (2016). Anticancer Res 36, 1,211–1,219.



---

**Generation and characterization of a MYF5 reporter human iPS cell line using CRISPR/Cas9 mediated homologous recombination.**

Wu J, Hunt SD, Xue H, Liu Y, Darabi R (2016). Sci Rep 6, 18759.



---

**Genetic modification in human pluripotent stem cells by homologous recombination and CRISPR/Cas9 system.**

Xue H, Wu J, Li S, Rao MS, Liu Y (2016). Methods Mol Biol 1307, 173–190.



---

**Generation and validation of PAX7 reporter lines from human iPS cells using CRISPR/Cas9 technology.**

Wu J, Hunt SD, Xue H, Liu Y, Darabi R (2016). Stem Cell Res 16, 220–228.



---

**A conserved C-terminal element in the yeast Doa10 and human MARCH6 ubiquitin ligases required for selective substrate degradation.**

Zattas D, Berk JM, Kreft SG, Hochstrasser M (2016). J Biol Chem 291, 12,105–12,118.



---

**Generation of human embryonic stem cell line expressing zsGreen in cholinergic neurons using CRISPR/Cas9 system.**

Zhou J, Wang C, Zhang K, Wang Y, Gong X, Wang Y, Li S, Luo Y (2016). Neurochem Res [published online ahead of print April 25, 2016]. Accessed June 14, 2016.



---

## 2015

---

### **Cloning-free CRISPR.**

Arbab M, Srinivasan S, Hashimoto T, Geijsen N, Sherwood RI (2015). *Stem Cell Reports* 5, 908–917.



---

### **Regulation of gene editing activity directed by single-stranded oligonucleotides and CRISPR/Cas9 systems.**

Bialk P, Rivera-Torres N, Strouse B, Kmiec EB (2015). *PLoS One* 10, e0129308.



---

### **Engineering human stem cell lines with inducible gene knockout using CRISPR/Cas9.**

Chen Y, Cao J, Xiong M, Petersen AJ, Dong Y, Tao Y, Huang CT, Du Z, Zhang SC (2015). *Cell Stem Cell* 17, 233–244.



---

### **CRISPR/Cas9-induced disruption of paraflagellar rod protein 1 and 2 genes in *Trypanosoma cruzi* reveals their role in flagellar attachment.**

Lander N, Li ZH, Niyogi S, Docampo R (2015). *MBio* 6, e01012



---

### **Metabolic engineering of *Escherichia coli* using CRISPR-Cas9 mediated genome editing.**

Li Y, Lin Z, Huang C, Zhang Y, Wang Z, Tang YJ, Chen T, Zhao X (2015). *Metab Eng* 31, 13–21.



---

### **Targeted disruption of *DNMT1*, *DNMT3A* and *DNMT3B* in human embryonic stem cells.**

Liao J, Karnik R, Gu H, Ziller MJ, Clement K, Tsankov AM, Akopian V, Gifford CA, Donaghey J, Galonska C, Pop R, Reyon D, Tsai SQ, Mallard W, Joung JK, Rinn JL, Gnirke A, Meissner A (2015). *Nat Genet* 47, 469–478.



---

### **Multiplex CRISPR/Cas9-based genome editing for correction of dystrophin mutations that cause Duchenne muscular dystrophy.**

Ousterout DG, Kabadi AM, Thakore PI, Majoros WH, Reddy TE, Gersbach CA (2015). *Nat Commun* 6, 6244.



---

### **Caspase-4 mediates non-canonical activation of the NLRP3 inflammasome in human myeloid cells.**

Schmid-Burgk JL, Gaidt MM, Schmidt T, Ebert TS, Bartok E, Hornung V (2015). *Eur J Immunol* 45, 2,911–2,917.





**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** 66 662 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 020 8328 2000

