

# Efficient editing of human iPS cells using vesicle-mediated delivery of Cas9-sgRNA protein complexes



Montse Morell<sup>1</sup>, Tatiana Garachtenko<sup>1</sup>, Lily Lee<sup>1</sup>, Mei Fong<sup>1</sup>, Thomas P Quinn<sup>1</sup>, Michael Haugwitz<sup>1</sup>, and Andrew Farmer<sup>1</sup> <sup>1</sup>Takara Bio USA, Inc., Mountain View, California, USA

## Abstract

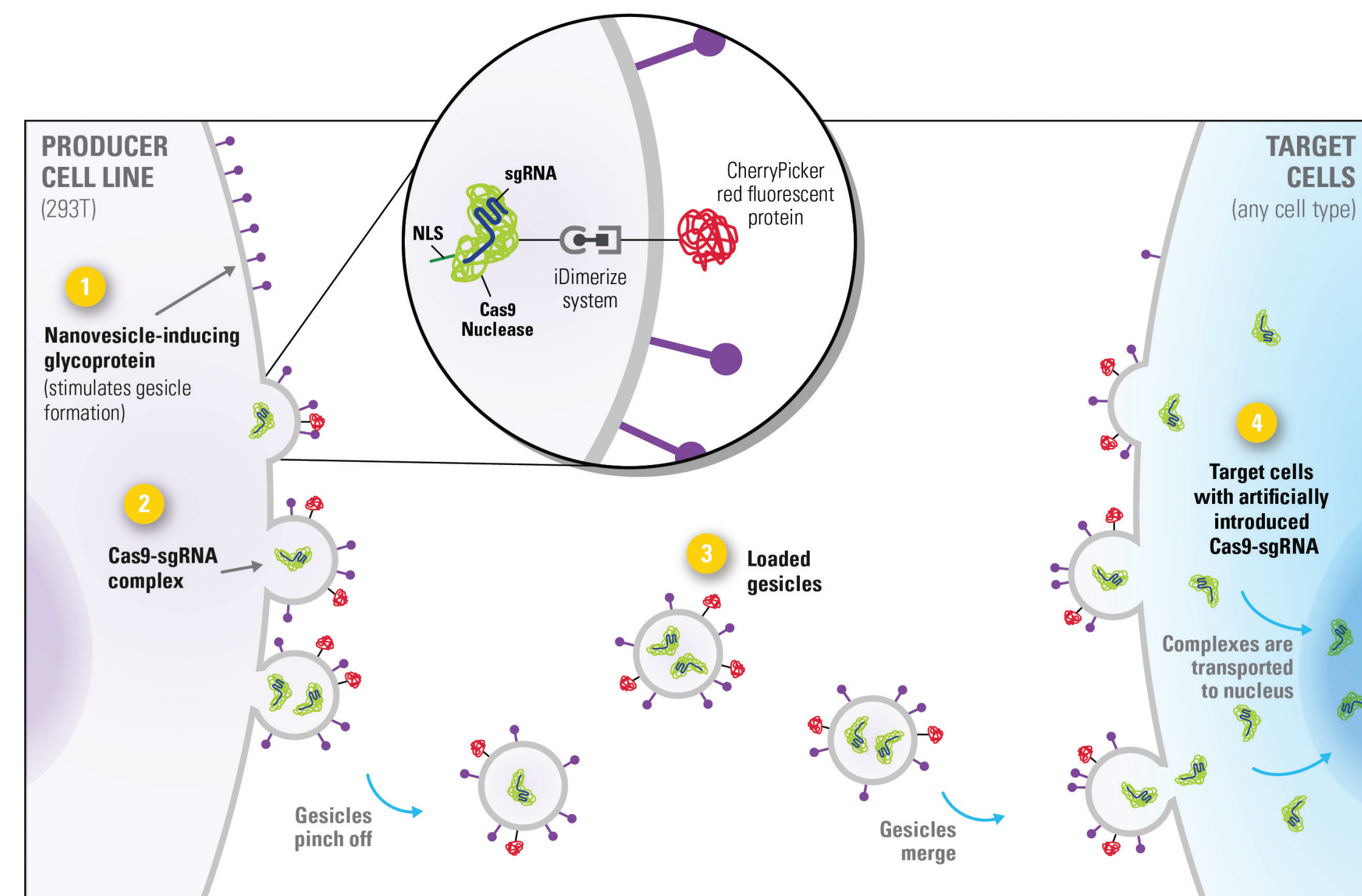
CRISPR/Cas9-based gene editing has revolutionized the field of cell biology. However, two significant challenges remain: obtaining efficient delivery of Cas9 to all cell types and achieving fewer off-target effects. It has been previously demonstrated that genome editing via direct delivery of Cas9-sgRNA ribonucleoproteins (RNPs) has the added benefit of decreased off-target effects due to the short duration of the RNP in the cell.

Here we report on a new RNP delivery method for footprint-free genome editing in a broad range of cell types, including human induced pluripotent stem cells (hiPS cells). Cas9 containing a nuclear localization signal (NLS) and an sgRNA of interest are simultaneously expressed in a mammalian packaging cell and actively packaged as an RNP complex into nanoparticles that originate at the plasma membrane. These cell-derived nanoparticles are called "vesicles". Vesicles are produced by the packaging cell due to the expression of a nanoparticle-inducing glycoprotein. We have adapted vesicle technology for the packaging of Cas9-sgRNA RNPs via a ligand-dependent dimerization system (iDimerize™ technology). This method enriches RNPs at the plasma membrane where they are incorporated into newly forming vesicles that are later harvested from the media. When added to target cells, vesicles will deliver the active RNP complexes into the cells.

In this work, we were able to demonstrate that vesicles carrying RNPs can mediate target-specific gene knockout in a broad range of cell types, including hiPS cells, without the toxic effects of other methods. In addition, this nanoparticle-based method allows for tight control of the timing and dose of the Cas9-sgRNA complexes, which decreases off-target effects. Due to the need to do genome editing in hiPS cells for the study of human biology and disease, we have optimized the use of vesicles to knock out genes in hiPS cells without affecting pluripotency.

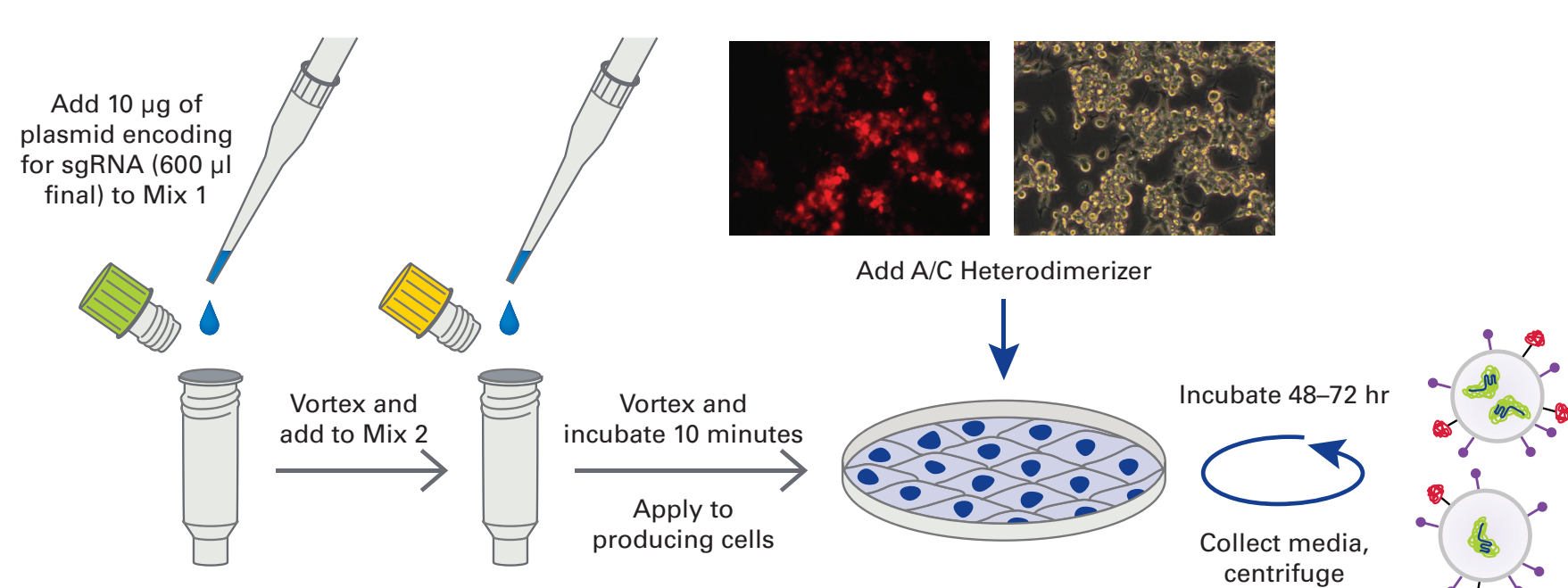
Overall, vesicles can be considered of high interest for genome editing, providing a direct, rapid, and transient method for delivering highly active Cas9-sgRNA RNPs into target cells.

## Mechanism of vesicle production for delivery of Cas9-sgRNA protein complex



**Production of nanovesicles (vesicles) for delivery of Cas9 protein together with sgRNA against a gene of interest.** Overview of Cas9 Vesicle production (Cat. #632613). (Step 1) In this illustration, expression constructs for Cas9 endonuclease, sgRNA against a target gene, CherryPicker™ red fluorescent protein, and other elements needed for vesicle production are cotransfected into the HEK 293T-based packaging cell line (Vesicle Producer 293T Cell Line, Cat. #632617). (Step 2) Overexpression of these proteins in the presence of the iDimerize ligand leads to the formation of vesicles containing sgRNA-loaded Cas9 protein, which are then collected from the media (Step 3). Addition of these vesicles to a target cell results in delivery of the Cas9-sgRNA complex to the nucleus and transient labeling of the membrane with CherryPicker protein (Step 4).

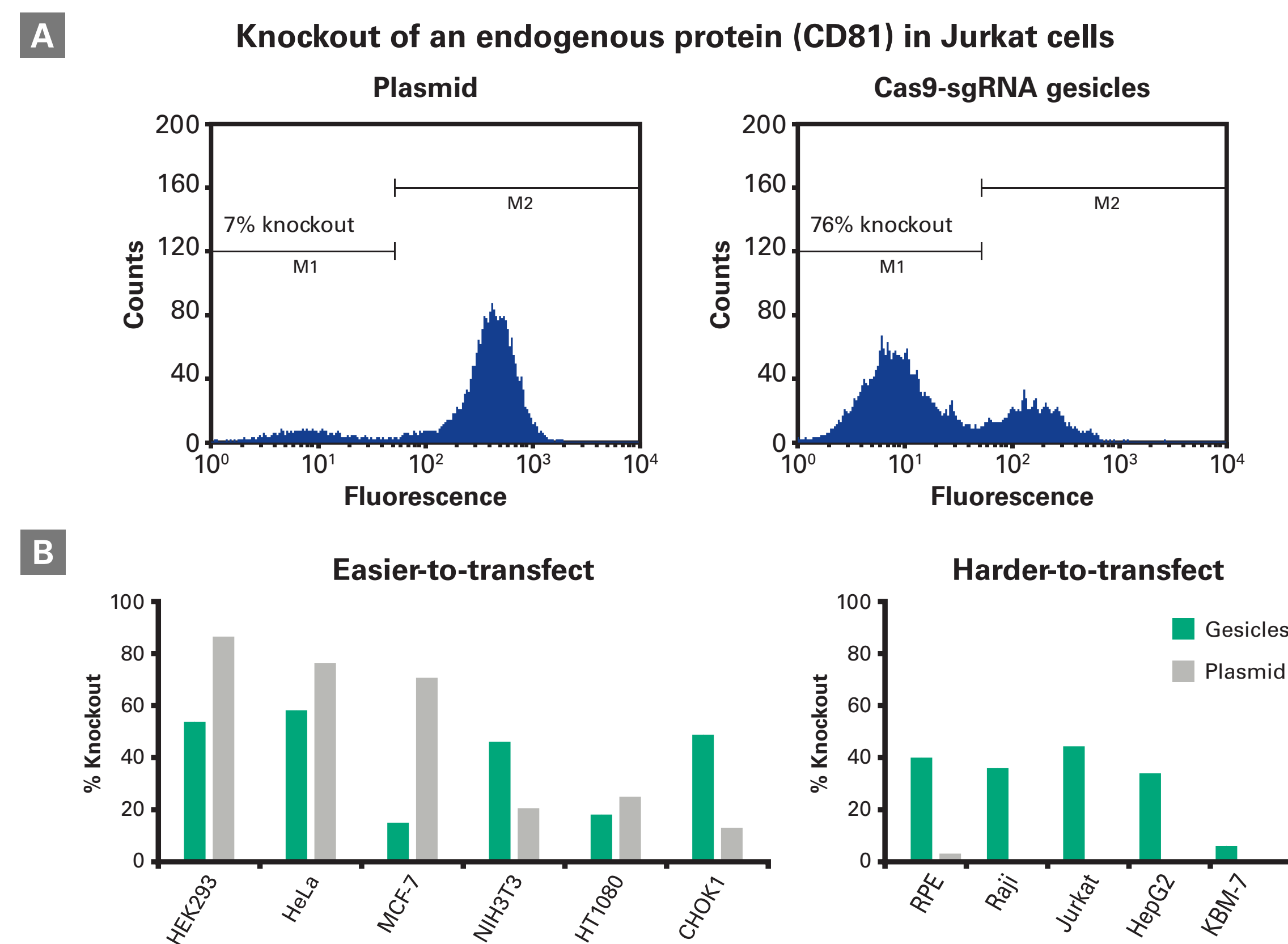
## Figure 1: Production of Cas9-sgRNA vesicles using the Guide-it™ packaging set



**Production of Cas9-sgRNA vesicles using the Guide-it packaging set.** The Guide-it vesicle packaging mix (Cat. #632616) contains lyophilized Xfect™ transfection reagent premixed with an optimized formulation of plasmids encoding for all the elements needed for vesicle production. Simply add your pGuide-it-sgRNA designed against your gene of interest, and then apply to Vesicle Producer 293T cells together with the iDimerize ligand. After 48-72 hr, centrifuge the cell media overnight at 7000 rpm. Decant the supernatant and resuspend the pellet (vesicles containing Cas9-sgRNA complexes). These Cas9 Vesicles are then ready to be added to target cells.

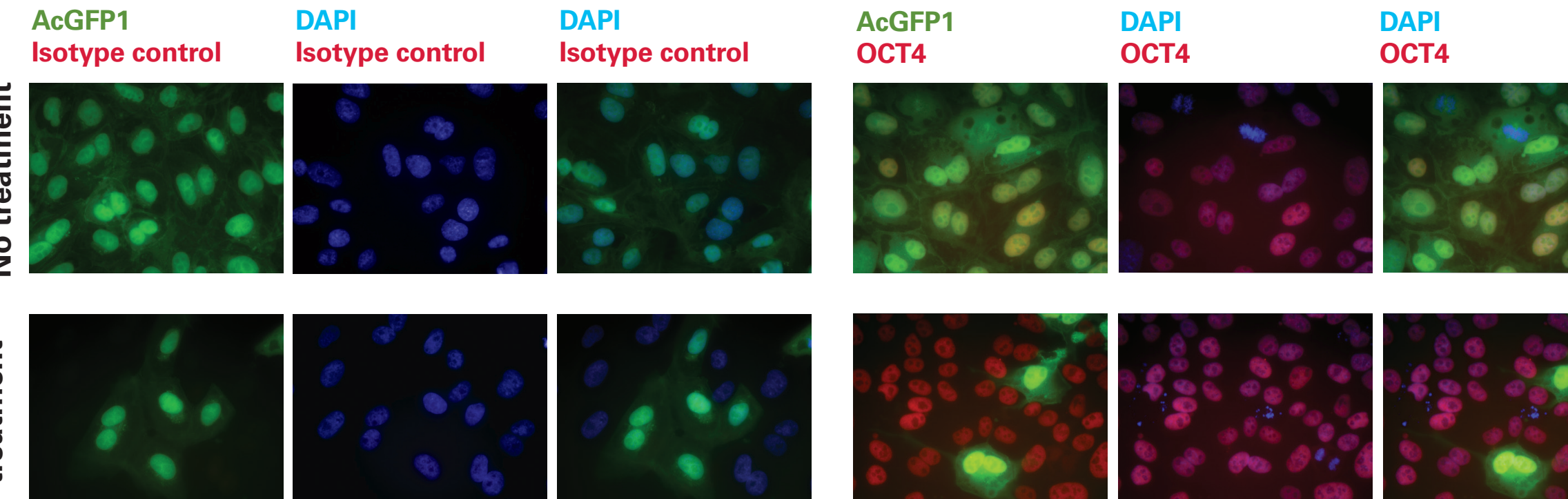
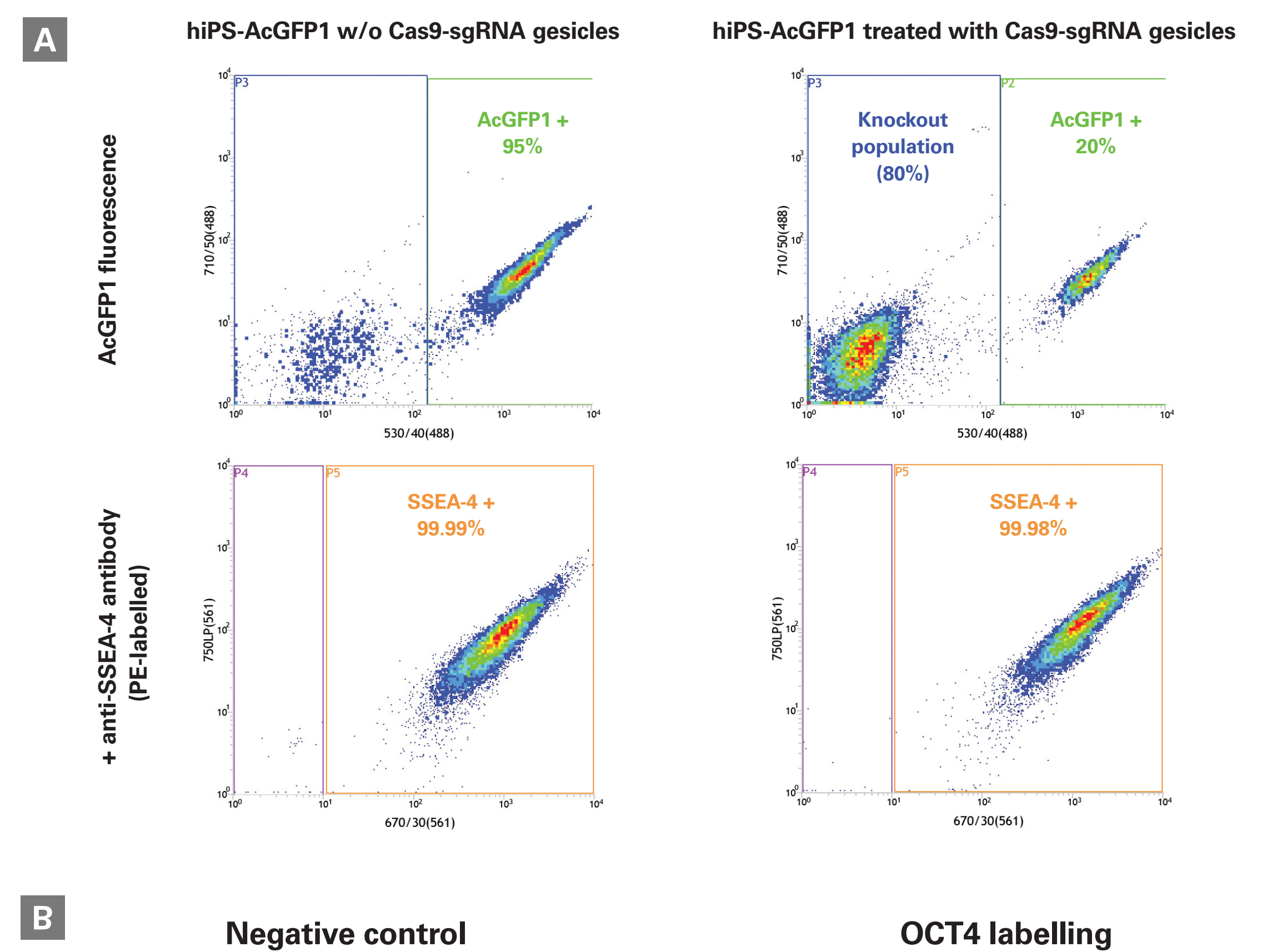
Clontech Laboratories, Inc. • A Takara Bio Company  
United States/Canada: +1 800.662.2566 • Asia Pacific: +1 858.919.7200 • Europe: +33 (0)1 3904 6880 • Japan: +81 (0)77 543 7267  
For Research Use Only. Not for use in diagnostic or therapeutic procedures. Not for resale. Takara® and the Takara logo are trademarks of Takara Holdings, Inc. Cellartis®, DEF-CS, and DEF-iPSC are trademarks of Takara Bio Europe AB. Clontech®, the Clontech logo, CherryPicker®, Guide-it®, iDimerize®, and Xfect are trademarks of Clontech Laboratories, Inc. All other trademarks are the property of their respective owners. Certain trademarks may not be registered in all jurisdictions. ©2016 Clontech Laboratories, Inc.

## Figure 2: Use of Cas9-sgRNA vesicles as a genome editing tool



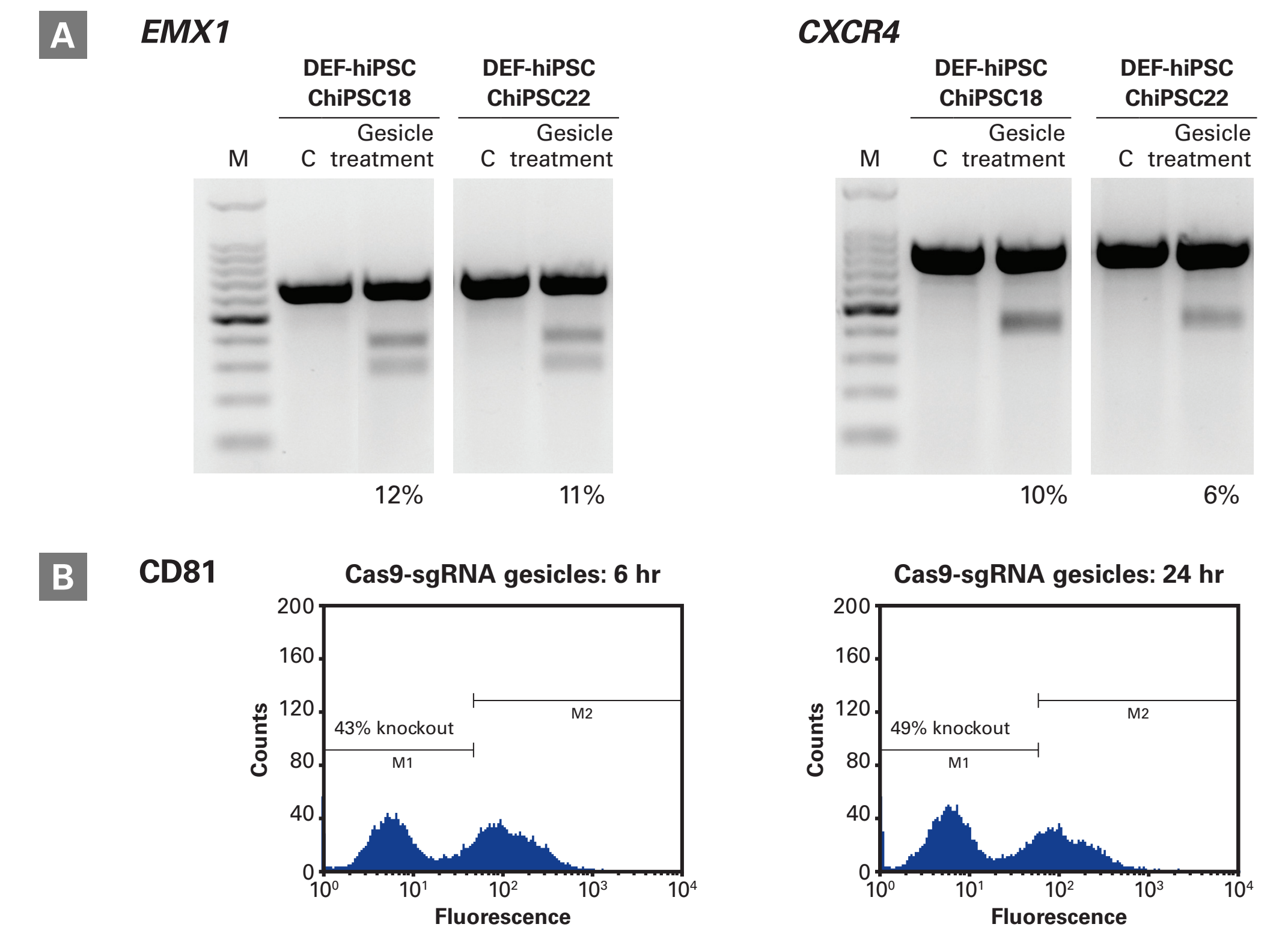
**Use of Cas9-sgRNA vesicles as a genome editing tool to knock out genes in multiple cell lines.** Panel A. The cell-surface protein receptor CD81 was knocked out in Jurkat cells using either plasmid cotransfection of Cas9 DNA and sgRNA, or Cas9-sgRNA vesicles (with the sgRNA against CD81 in both instances). Knockout efficiencies were measured six days later via antibody labeling of CD81 followed by flow cytometry. The knockout efficiency obtained via vesicles was significantly higher than the results achieved with plasmid transfection. Panel B. Cas9-sgRNA vesicles outperform plasmid transfection in harder-to-transfect cell types. Cell lines were created that contained an integrated ZsGreen1 fluorescent protein expression cassette. In this system, gene editing of ZsGreen1 can be monitored by a loss of green fluorescence. These cell lines were treated with Cas9-sgRNA vesicles (with the sgRNA against ZsGreen1), and then analyzed by flow cytometry. Cas9-sgRNA protein complex delivery and ZsGreen1 knockout via vesicles was efficient and comparable to plasmid-based delivery in easier-to-transfect cell types (left graph) and surpassed the results achieved via plasmid-based delivery in harder-to-transfect cell types (right graph).

## Figure 3: Knockout of a fluorescent protein in hiPS cells



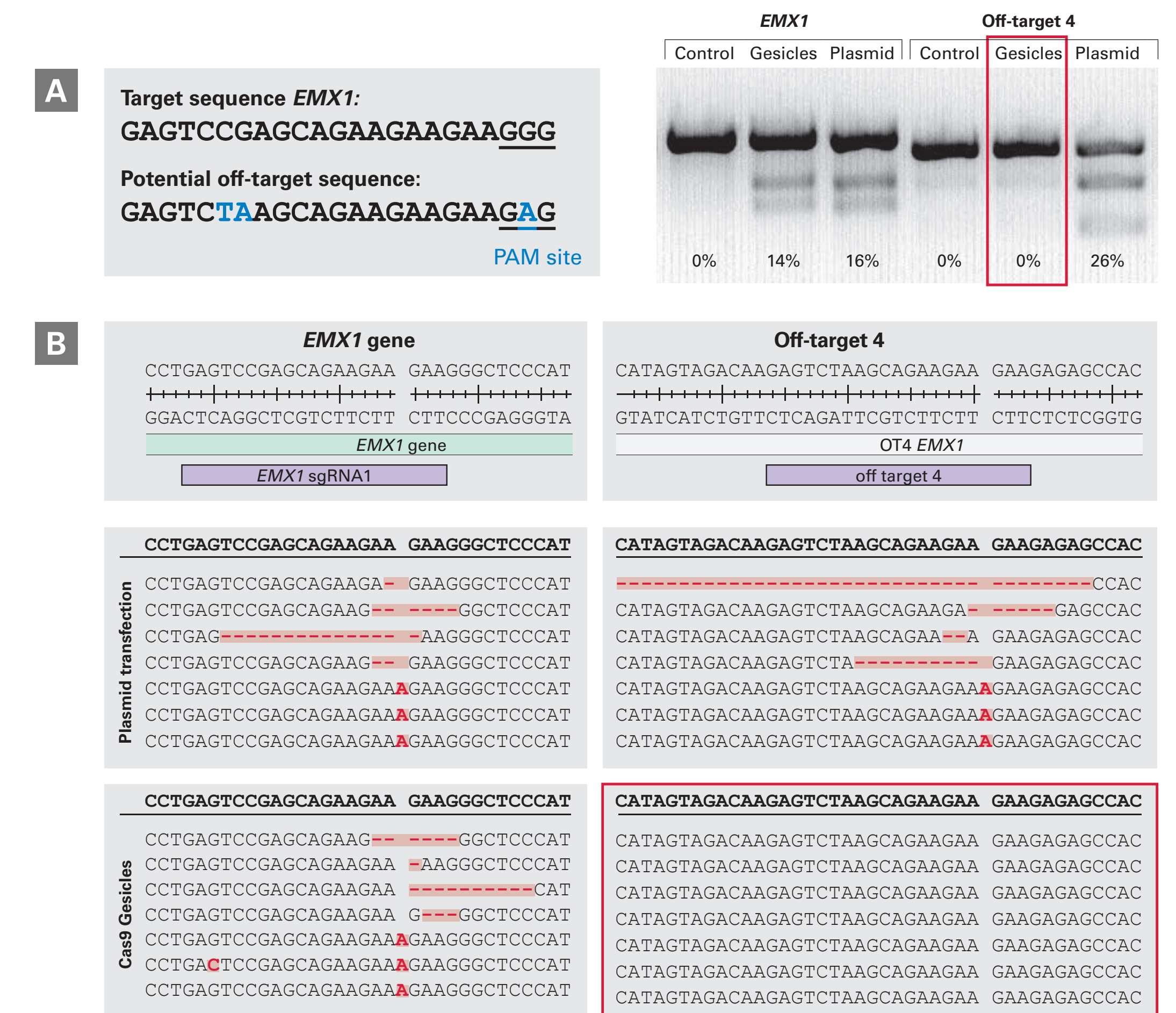
**Knockout of a fluorescent protein in hiPS cells.** Panel A. Cas9-sgRNA vesicles targeting AcGFP1 were added to Cellartis® Human iPS Cell Line 22 (DEF-iPSC™ ChiPSC22, Cat. #Y00325) stably expressing AcGFP1, cultured under non-differentiating conditions using the Cellartis DEF-CS™ Culture System (Cat. #Y30010). After vesicle treatment, AcGFP1 expression was determined via flow cytometry in untreated DEF-iPSC ChiPSC22-AcGFP1 (top left) and vesicle-treated DEF-iPSC ChiPSC22-AcGFP1 (top right). Both samples were also labelled with anti-SSEA4 (PE) antibody (bottom). Vesicle-mediated delivery of Cas9-sgRNA complexes induced knockout in about 75% of cells without a loss of pluripotency. Panel B. Immunohistochemistry was performed on the DEF-iPSC ChiPSC22-AcGFP1 cells six days after vesicle addition. Cells were permeabilized and stained with an antibody against OCT4 (eFluor 570-labelled) or an IgG2a isotype control eFluor 570. Cells treated with the vesicles presented a mixed population of fluorescent and non-fluorescent cells (a product of successful gene editing of the AcGFP1 gene by Cas9 Vesicles), all of them positive for OCT4 expression. Therefore, vesicle treatment left the pluripotency of the edited cells unaffected.

## Figure 4: Knockout of endogenous genes in hiPS cells



**Efficient knockout of endogenous genes in hiPS cells using vesicles.** Panel A. The genes EMX1 (left) and CXCR4 (right) were knocked out using Cas9-sgRNA vesicles in DEF-iPSC ChiPSC18 and DEF-iPSC ChiPSC22 cell lines. Knockout efficiencies were determined using a Guide-it Resolvase assay (Guide-it Mutation Detection Kit, Cat. #631443). A control sample that was not treated with Cas9 was included for comparison (Control, C). The percentage of DNA cleavage was determined by densitometry. The knockout could be detected in all the samples. Panel B. Successful knockout of CD81 in hiPS cells. Vesicles containing Cas9-sgRNA complexes designed to target human CD81 were harvested and added to Cellartis Human iPS Cell Line 18 (Cat. #Y00305), cultured with the DEF-CS Culture System for six or 24 hr, and then cultured in vesicle-free DEF-CS culture media for an additional seven days. The surface expression of CD81 on vesicle-treated cells was determined via flow cytometry. Graphs show DEF-iPSC ChiPSC18 cells after six hr (left) or 24 hr (right) of vesicle treatment, labeled with anti-CD81 (FITC) antibodies. There is an increase in knockout percentage when the incubation time is increased.

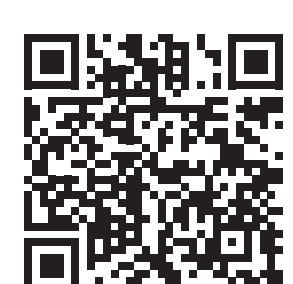
## Figure 5: Cas9 vesicles reduce off-target effects



**The use of Cas9 Vesicles reduces off-target effects compared with plasmid-based treatment.** Panel A. Decrease of off-target effects detected via Guide-it Resolvase assay. HEK 293T cells were either simultaneously transfected with plasmids encoding Cas9 and a sgRNA against EMX1, or treated with Cas9-sgRNA vesicles. After 72 hr, the EMX1 gene and a potential off-target locus (off-target 4) were amplified and the percentage of DNA cleavage was determined using a Guide-it Resolvase assay. With the Cas9 Vesicles, no off-target effect could be detected. Panel B. Indels generated after Cas9 targeting of EMX1. Indels were identified using the Guide-it Indel Identification Kit (Cat. #631444). Sequencing data for the different clones revealed a range of deletions and insertions (highlighted in red) in both sites (target site EMX1 as well as off-target 4) in the case of cotransfection with Cas9 and sgRNA plasmids. In the case of the cells treated with Cas9 Vesicles, indels could only be detected at the target site; the off-target site was not mutated.

## Conclusions

- Vesicles provide a highly efficient means of producing, packaging, and directly delivering Cas9 protein complexed with sgRNA to any target cell, including hiPS cells
- Delivery of Cas9 via vesicles leads to levels of genome modification similar to plasmid delivery, but with the added benefit of drastically reduced editing of off-target sites
- Vesicles can be used to knock out genes in hiPS cells without any change in cell pluripotency



Scan to download your copy of this poster and view additional posters, or visit <http://info.clontech.com/CSHL-2016-Poster-Signup.html>

800.662.2566

Visit us at [www.clontech.com/vesicle](http://www.clontech.com/vesicle)

Clontech Takara cellartis