

Product Information

Human RPTEC OCT2 KO Cells

Catalog Number **MTOX1080**

Storage Temperature $-130\text{ }^{\circ}\text{C}$ or below
in liquid nitrogen vapor phase

TECHNICAL BULLETIN

Product Description

CompoZr® zinc finger nuclease (ZFN) technology is a fast and reliable way to manipulate the genome in a targeted fashion. ZFNs are naturally occurring proteins that can be engineered to bind DNA at a sequence-specific location and create a double strand break (www.sigma.com/zfn). The cell's natural machinery repairs the break in one of two ways: non-homologous end joining or homologous recombination. The non-homologous end joining pathway typically produces small modifications (indels) at the targeted locus that may result in a functional knockout. Single cell clones are then isolated, tested for the desired modification, and expanded to establish stable cell lines.

Kidney toxicity is a major concern during drug development. Although primary human kidney cells are available and a handful of immortalized kidney cell lines currently exist, there are concerns with passage limitations and limited functionality in these systems. Human primary proximal tubule epithelial cells (RPTEC) were modified with ZFNs to effectively extend cell proliferation. The resulting OCT 2 Knockout (KO) cell line was characterized for the presence of proximal tubule cell markers as well as several functional properties, including response to several known human nephrotoxics similar to human primary cells and robust expression, and activity of key renal uptake and efflux transporters including OCT2, OATP4C1, OAT1, OCTN1, MRP2, MRP4, P-gp, MATE1, MATE2-K, PEPT1, and PEPT2.

This product consists of ZFN engineered RPTEC OCT2 KO Cells. They are intended for use with SA7K Clone Control Cells (Catalog Number MTOX1030) for a wide variety of kidney cell based assays.

Components

Human RPTEC OCT2 KO Cells 1 vial
Catalog Number MTOX1080
Vial contains $\geq 3 \times 10^6$ modified RPTEC cells.

Note: Neither media nor supplements are supplied with the vials. These must be obtained prior to receiving the vials.

The cryoprotectant medium used is Cell Freezing Medium-DMSO 1 \times , Catalog No. C6164

Parental Cell Line: Zen Bio HK016

Note: Please see product datasheet from Zen Bio for additional information about the origin of these cell lines. Cytogenetic information is based on initial seed stock at Sigma Life Science. Cytogenetic instability has been reported in the literature for some cell lines.

Cell Line Description

Clone: RPTC-HK016 Zen Bio
Organism: *Homo sapiens* (human)

Tissue: Renal Proximal Convoluted Tubule Epithelial Cells

Age: 8.8 years

Gender: Female

Ethnicity: African American

Morphology: Epithelial

Growth properties: Adherent

Species-specific PCR Evaluation:

The cells were confirmed to be of human origin and no mammalian interspecies contamination was detected.

PCR Evaluation for *Mycoplasma* species contamination:
Negative

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

Storage/Stability

Upon receiving a shipment of frozen cells it is important the end user gives the shipment attention without delay. To ensure the highest level of viability, thaw the vial and initiate the culture as soon as possible upon receipt. If upon arrival, continued storage of the frozen culture is necessary, it should be stored in liquid nitrogen vapor phase and not at $-70\text{ }^{\circ}\text{C}$. Storage at $-70\text{ }^{\circ}\text{C}$ will result in loss of viability.

Procedures

Protocol for Thawing and Seeding 24 Well Plates

Note: One cryovial of Human RPTEC OCT2 KO Cells contains enough cells to seed one plate.

Reagents and Equipment Required but Not Provided for Thawing and Seeding

Note: Neither media nor supplements are supplied with the vials. These must be obtained prior to receiving the vials.

- RPTEC Complete Supplement (Catalog Number MTOXRCSUP – 30 mL)
- RPTEC Tox Supplement (Catalog Number MTOXRTSUP – 6 mL)
- Minimum Essential Medium Eagle Alpha Modification Medium ($\text{MEM}\alpha$) (Catalog Number M4526)
- L-glutamine (Catalog Number G7513)
- Gentamicin solution (Catalog Number G1397)
- Amphotericin B solution (Catalog Number A2942)
- HBSS (Catalog Number H6648)
- BSL-2 hood
- Cell culture incubator

Medium Preparation

Prepare Complete Medium by supplementing 500 mL of $\text{MEM}\alpha$ (Catalog Number M4526) with 6.25 mL of L-glutamine (Catalog Number G7513), 30 mL of RPTEC Complete Supplement (Catalog Number MTOXRCSUP), 0.3 mL of Gentamicin solution (Catalog Number G1397), and 0.03 mL of Amphotericin B solution (Catalog Number A2942). Complete Medium can be stored at $2\text{--}8\text{ }^{\circ}\text{C}$ for up to 1 month. This medium is formulated for use with a 5% CO_2 in air atmosphere.

Prepare Tox Medium by supplementing 500 mL of $\text{MEM}\alpha$ (Catalog Number M4526) with 6.25 mL of L-glutamine (Catalog Number G7513), 6 mL of RPTEC Tox Supplement (Catalog Number MTOXRTSUP), 0.3 mL of Gentamicin solution (Catalog Number G1397), and 0.03 mL of Amphotericin B solution (Catalog Number A2942). Tox Medium can be stored at $2\text{--}8\text{ }^{\circ}\text{C}$ for up to 1 month. This medium is formulated for use with a 5% CO_2 in air atmosphere.

Thawing of Frozen Cells

1. Thaw the vial by gentle agitation in a $37\text{ }^{\circ}\text{C}$ water bath. To reduce the possibility of contamination, keep the O-ring and cap out of the water. Thawing should be rapid (~ 2 minutes).
2. Remove the vial from the water bath as soon as the contents are thawed, and decontaminate by dipping in or spraying with 70% ethanol. All of the operations from this point on should be carried out under strict aseptic conditions.
3. Transfer the vial contents to a centrifuge tube containing 9.0 mL of RPTEC Complete Medium and spin at $125 \times g$ for 3–5 minutes.
4. Resuspend cell pellet with the RPTEC Complete Medium. It is important to avoid excessive alkalinity of the medium during recovery of the cells. It is suggested, prior to the addition of the vial contents, the culture vessel containing the REPTC Complete Medium be placed into the incubator for at least 15 minutes to allow the medium to reach its normal pH (7.0–7.6) and temperature ($37\text{ }^{\circ}\text{C}$).

Transporter Assays

1. Resuspend cell pellet in 12 mL of RPTEC Complete Medium.
2. Add 0.5 mL of suspension per well of 24 well plate. Cells should be greater than 95% confluent after 2 days.

Uptake/Efflux Transporter Assays

1. For assays involving the use of chemical inhibitors, add the desired concentration of the inhibitor to selected wells and pre-incubate for 30 minutes.
2. Following the pre-incubation with inhibitors, add desired concentration of substrates and incubate for 30 minutes.
3. Stop the reaction by washing the cells three times with ice-cold HBSS buffer.
4. Prepare and analyze the samples according to established protocols.

Toxicity Assay

1. Resuspend cell pellet in 12 mL of RPTEC Complete Medium and add 0.5 mL per well of 24 well plate. Cells should be greater than 95% confluent after 2 days.
2. On day of treatment, change medium to Tox Medium. Prepare and dilute compound to generate a dose concentration response curve. Treat cells with appropriate volume of compound to obtain final concentrations.
3. After 48–72 hours, determine toxicity of compound. Note: The use of MTT Based *In Vitro* Toxicology Assay Kit (Catalog Number TOX1) is recommended.

CompoZr is a registered trademark of Sigma-Aldrich Co. LLC.

ANA,TL,MAM 01/18-1

These products are covered by the “Consumable Vial” License as described in Exhibit 1.

**EXHIBIT
1
LICENSE AGREEMENT - ADME/TOX CELL
LINES**

This Product and its use are the subject of one or more of the following patents controlled by Sangamo BioSciences, Inc.: U.S. Patent Nos. 6,534,261, 6,607,882, 6,746,838, 6,794,136, 6,824,978, 6,866,997, 6,933,113, 6,979,539, 7,013,219, 7,030,215, 7,220,719, 7,241,573, 7,241,574, 7,585,849, 7,595,376, 6,903,185, 6,479,626, US20030232410, US20090203140 and corresponding foreign patent applications and patents.

BEFORE OPENING OR USING THIS PRODUCT, PLEASE READ THE TERMS AND CONDITIONS SET FORTH IN THIS LICENSE AGREEMENT. YOUR USE OF THIS PRODUCT SHALL CONSTITUTE ACKNOWLEDGMENT AND ACCEPTANCE OF THESE TERMS AND CONDITIONS. If you do not agree to use this Product pursuant to the terms and conditions set out in this License Agreement, please contact Sigma Technical Services within ten days of receipt to return the unused and unopened Product for a full refund; provided, however, that custom-made Products may not be returned for a refund.

The purchase of this Product conveys to you, the buyer, the non-transferable right to use the purchased Product for Licensed Research Use (see definition below) subject to the conditions set out in this License Agreement. If you wish to use this Product for any purpose other than Licensed Research Use, you must first obtain an appropriate license (see information set out below)

This Product may not be used for any purpose other than Licensed Research Use. Product, as used herein, means the Cell Line and any derivative cells or cell lines created by the buyer which contain and/or incorporate genetic information derived from the Cell Line. Your right to use this Product for Licensed Research Use is subject to the following conditions and restrictions:

The type of license you have subject to this Agreement (Evaluation, Annual, Extended and Consumable) is listed on the outside of the package and/or on the invoice you received from us.

1. “Licensed Research Use” means any use for research purposes, other than:

- (a) Licensing, selling, distributing, or otherwise providing the product or modified versions of it to any third party other than Sigma and its affiliates as provided herein: provided however, that you may provide the product or modified versions of it to researchers within your research organization located at the same research facility or campus who are similarly bound to the use restrictions herein;
- (b) GMP production of therapeutic, diagnostic, prophylactic or other medicinal products intended for use in humans or non-human animals, or any other industrial use solely to the extent involving commercial sale of a product or service. If a molecule or any

derivative of such molecule is used in or administered to humans, then the production of such molecule shall be deemed to be GMP production and therefore in violation of this License Agreement;

2. You may not transfer the Product, its components, or any materials made through the use of this Product, including further modified cells, to any third party without prior written approval of Sigma. Notwithstanding the foregoing, the Product or materials made through use of the Product may be transferred by you without such prior written approval to your legal affiliates or bona fide third party contractors performing paid work on your behalf, provided the use by such third party contractors is limited to performance of work for you and any results or product of such work shall not be shared by the third party contractor with any other person.

3. For Products covered by an "Evaluation" License, this Agreement shall expire ninety (90) days following your receipt of the Product.

4. For Products covered by an "Annual" License, this Agreement shall expire three hundred and sixty-five (365) days following your receipt of the Product.

5. For Products covered by an "Extended" License, this Agreement shall remain in force in perpetuity following your receipt of the Product.

6. For Products covered by a "Consumable" License, you may:

(a) Not expand, propagate, seed, bank, cryopreserve or store the product, its components, or any materials made through the use of this Product, outside of the original shipping device, including but not limited to vial, transwell plate and Petaka cell culture device. The Product can only be transferred and/or incorporated in an end-use assay. All Products that are not utilized in an end-use assay shall be immediately destroyed.

(b) Use Products for the provision of commercial services to a third party for monetary gain, but limited to only delivery of research results and data from use of the Product, and not the Product, its components, or any materials made through the use of this Product, including further modified cells.

7. Products covered by a "Consumable Vial" License can only be incorporated in an end-use assay, seeded directly on the final assay plate (including, but not limited to 6-well, 12-well, 24-well, 48-well, or 96-well plates). All Products that are not utilized in an end-use assay plates shall be immediately destroyed.

8. Products delivered in a Petaka cell culture device can only be incorporated in an end-use assay for up to two (2) transwell plates (including, but not limited to 24-well or 96-well plates). All Products that are not utilized in an end-use assay for up to two (2) transwell plates shall be immediately destroyed.

9. Your right to use the Product will terminate immediately upon expiration of this License Agreement, or in the event that you fail to comply with these terms and conditions. You shall, upon such termination of your rights, destroy all Product, any modified versions of the Product, and components thereof in your control, and notify Sigma of such in writing.

For information on purchasing a license to this Product for purposes other than Licensed Research Use, contact your local Sigma Sales representative, who will refer you to the proper licensing representative, or in the USA call 800-325-3010.