

TransfeX[™] Transfection of Plasmid DNA into Primary Human Mammary Epithelial Cells (HMECs)

HMECs (ATCC[®] Cat. No. PCS-600-010[™]) are primary, normal, human, epithelial cells isolated from adult mammary gland tissue. ATCC has achieved transfection efficiencies of approximately **60%** using the protocol described below.

General Considerations for using the TransfeX transfection reagent:

- All steps should be performed in a biosafety cabinet using proper aseptic technique.
- **Cell conditions.** Cells should be passaged at least once after thaw and the use of lowpassage cells is recommended. Passage the cells 18-24 hours before transfection to ensure the cells are actively dividing and that they will be at the appropriate cell density at the time of transfection. Make sure that the cells are healthy and are ≥ 90% viable, prior to transfection.
- Seeding density. Cell density should be 50-80% confluent on the day of transfection. See specified seeding density in the individual protocols and in Table 1. *Note: Determine the optimal cell density for each cell type in order to maximize transfection efficiency.*
- **DNA purity.** Use highly purified plasmid preps that are free from phenol or other contaminants. Plasmid DNA preps that are endotoxin-free are desirable.
- **Presence of antibiotics and other inhibitors.** Antibiotics will inhibit transfection complex formation and therefore should be excluded from the complex formation step. Transfection complexes can be added to cells grown in complete culture medium containing serum and low levels of antibiotics if required.
- **Complex formation conditions.** Prepare TransfeX Reagent and DNA complexes in serumfree growth medium. ATCC recommends using Opti-MEM I Reduced-Serum Medium to dilute the DNA before complex formation.

Materials required:

alog No.
C [®] PCS-600-010™
C [®] PCS-600-030
C [®] PCS-600-040
C [®] ACS-4005
Technologies™ 31985-062



Protocol:

The following protocol describes how to transfect plasmid DNA into primary HMECs using the TransfeX Reagent in **a single well of a 24 well plate.** The reaction may be scaled up at needed. Please refer to Table 1 for recommended reaction conditions for other dish or plate sizes.

A. Preparation of the cells for transfection

The day before transfection:

1. Count and measure cells for density and viability.

Plate **25,000** cells per well in complete growth medium (**Mammary Epithelial Cell Basal Medium supplemented with the Mammary Epithelial Cell Growth Kit**). Cell density should be **75%** confluent on the day of transfection.

2. Incubate cells overnight at 37°C with 5% CO₂.

The day of transfection:

- 1. Remove old media.
- 2. Replace old media with fresh complete growth media (excluding Heparin Sulfate) to a total volume of 0.5 mL/well.

B. Preparation of the DNA:TransfeX transfection complexes

- 1. Warm TransfeX, plasmid DNA, and Opti-MEM I Reduced-Serum Medium to room temperature and vortex gently to mix.
- 2. Pipette 50 µL Opti-MEM I Reduced-Serum Medium into a sterile microcentrifuge tube.
- 3. Add 0.5 μL (1.0 μg/μL) plasmid DNA.
- 4. Mix thoroughly with gently pipetting.
- 5. Add 1 µL TransfeX Reagent to the diluted DNA mixture. Note: Do not let the pipette tip or the reagent come into contact with the sides of the plastic tube.
- 6. Mix TransfeX:DNA complexes thoroughly using either a vortex or by pipetting briefly.
- 7. Collect contents at bottom of the tube using a mini-centrifuge.
- 8. Incubate TransfeX:DNA complexes at room temperature for 15 minutes.

C. Addition of DNA:TransfeX transfection complexes to the wells

- 1. Distribute the complexes to the cells by adding **50 μL** of the complexes drop-wise to different areas of the wells.
- 2. Gently rock the culture vessel back and forth and from side to side to evenly distribute the TransfeX:DNA complexes.

D. Post-Transfection Handling

- 1. Incubate for **24-72** hours. Replace transfection medium with fresh complete growth medium every 24 hours post transfection.
- 2. Wait for 18-24 hours post-transfection before assaying for transgene expression



Table 1: Recommended Reaction Conditions for different size culture vessels.

Culture Vessel	96 well plate	24 well plate	12 well plate	6 well plate	10 cm dish
Surface area	0.35 cm ²	1.9 cm ²	3.8 cm ²	9.6 cm ²	59 cm ²
Cell seeding	1.2 x 10 ⁴ per well	4.8 x 10 ⁴ per well	9.6 x 10 ⁴ per well	19 x 10 ⁴ per well	114 x 10 ⁴ per well
Complete Growth Medium	0.1 mL	0.5 mL	1.0 mL	2.0 mL	10 mL
Opti-MEM I Reduced Serum Medium	20 µL	50 µL	100 µL	200 µL	1.0 mL
DNA (1 µg/µL stock)	0.1 µg	0.5 µg	1.0 µg	2.5 µg	15 µg
TransfeX Reagent	0.2 µL	1.0 µL	2.0 µL	5.0 µL	30 µL

Notes:

Include a control condition consisting of an empty vector plasmid or a plasmid expressing GFP.