

Data Sheet • 293Ta Lentiviral Packaging Cell Line • Catalog No. CLv-PK-01

Description

293Ta packaging cells for lentivirus production using optimized packaging plasmids with lentiviral vector-based constructs.

Contents: 1 ml of 293Ta lentiviral packaging cells (1.5x10⁶ cells/ml). Low passage and authenticated. Cell contains: Adeno and SV-40 viral DNA sequences Growth properties: Adherent Organism: Homo sapiens (human) Source organ: Kidney Morphology: Epithelial

Growth medium: DMEM supplemented with 10% fetal bovine serum and Penicillin/Streptomycin Temperature: 37.0°C

Thawing and Plating 293Ta Cells

- 1. Remove the vial from liquid nitrogen freezer and immediately place it into a 37°C water bath. Remove the vial before the last bit of ice disappears to avoid overheating the cells. Wipe the vial with 70% ethanol.
- 2. Transfer thawed cells to a sterile 15ml or 50ml tube containing ~10 ml of complete growth medium (either cold or room temperature). Gently swirl the tube to mix.
- 3. Spin the cell suspension for 5 minutes at 300xg.
- Carefully discard supernatant as completely as possible. 4.
- 5. 6. Tap the tube gently to dislodge and loosen the cell pellet.
- Add ~10 ml of complete growth medium, and swirl the tube to suspend the cells.
- 7. Transfer the cell suspension to a 10cm culture dish or other appropriate culture vessel. The culture dish or vessel must be plasma tissue-culture treated.

Important Note: The 293Ta packaging cells were frozen in a DMSO containing medium. DMSO is toxic to cells and must be washed off before seeding the cells into a culture vessel. If DMSO is not removed, the cells will not attach and massive cell death will follow. If the culture dish is not plasma tissue-culture treated, the cells will not attach either.

Passaging

- 1. Discard culture medium.
- 2. Briefly rinse the cell layer with PBS.
- 3. Add 0.05% trypsin-EDTA solution to dish (just enough to cover the bottom of dish) and observe cells under an inverted microscope until the cell layer is dispersed.
- 4. Add >3x volume of complete growth medium (For example, add 9 ml or more of complete growth medium if 3 ml of trypsin-EDTA solution is used in step 3). Aspirate cells by gentle pipetting.
- 5. Transfer appropriate aliquots of the cell suspension to new culture vessels.
- 6. Incubate cells at 37°C with 5% CO₂.

Subcultivation ratio: Subcultivation ratio of 1:4 to 1:8 is recommended Medium renewal: Every 2 to 3 days Freeze medium: Complete growth medium supplemented with 8% (v/v) DMSO Storage temperature: Liquid nitrogen vapor phase

GeneCopoeia, Inc. 9620 Medical Center Drive, #101 Rockville, Marvland 20850 Tel: 301-762-0888 Fax: 301-762-8333 Email: inquiry@genecopoeia.com Web: www.genecopoeia.com

GeneCopoeia Products are for Research Use Only