

# Addressing Large-Scale Therapeutic Virus Production Using High Quality Grade PEI-Based Transfection Reagents

Mathieu Porte, Alengo Nyamay’Antu, Marine Ricordel, Mégane Denu, Patrick Erbacher

Polyplus-transfection\*, Vectura, 75 rue Marguerite Perey, 67400 Illkirch, France

## 1. Introduction

Gene and cell therapy-based medicines are experiencing resurgence due to the introduction of “next generation” transfer viral vectors, which have demonstrated improved safety and efficacy. Adeno Associated Virus (AAV) and Lentivirus are very commonly used in therapeutics and often produced using PEI-mediated transient transfection in HEK-293 or HEK-293T cells. The critical raw materials needed for cGMP vector production must be sourced from approved suppliers and should have gone through a rigorous testing program to reduce the risk of introducing adventitious agents into the production process. Polyplus-transfection now provides PEIpro®, the unique PEI-based transfection reagents available in different quality grades, allowing a seamless transition from process development with PEIpro®-HQ to cGMP biomanufacturing with PEIpro®-GMP. Here, we describe an optimized PEI-based virus production process for high-yielding viral vector production, compatible with different cell culture adherent and suspension systems. We further demonstrate the robust viral vector production yields, as well as the adaptability and reliability of the PEI-based transient gene expression approach to efficiently manufacture GMP-grade viral vectors at a sufficiently large scale for more advanced clinical trials, and in fine to drive commercialization of therapeutic vectors.

## 2. Optimized Transient Transfection for Virus Production

Figure 1: Optimization Process of PEI Polymer Chemistry

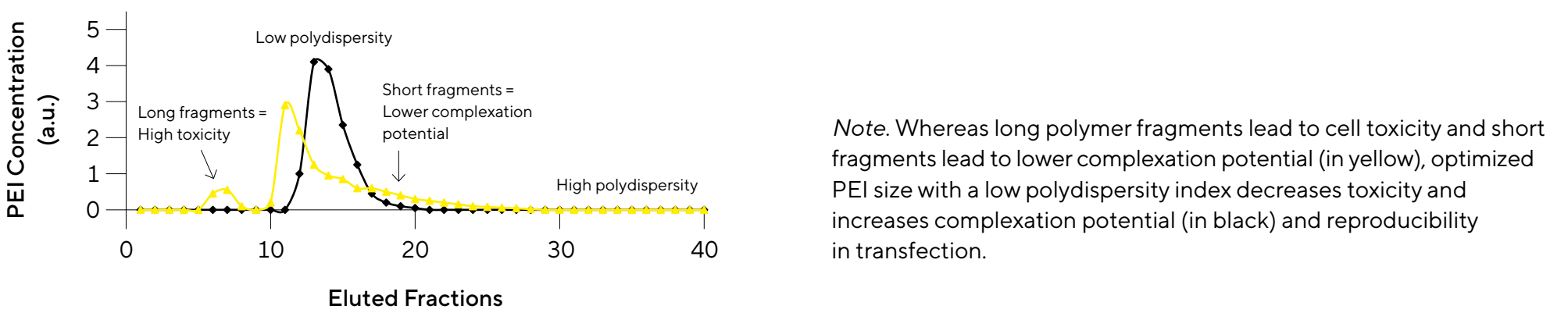
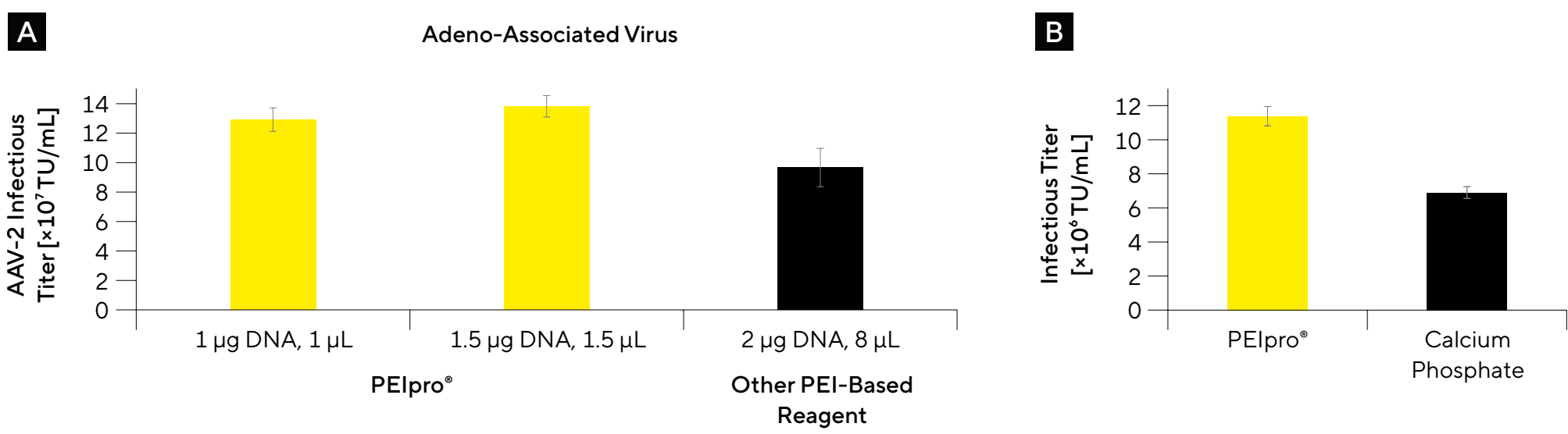


Figure 2: PEIpro® Produces More Virus With Less Reagent and Lower DNA Amount Compared to Another PEI-Based Reagent and Calcium Phosphate Transfection



Note. (A) Suspension HEK-293T cells were seeded at  $1 \times 10^6$  cells/mL in serum-free medium and transfected with PEIpro® and another PEI-based reagent following the recommended protocols. AAV-2 were produced with Helper Free Packaging System (Cell Biolabs, San Diego, CA) and titers were measured 72 hr after transfection using a GFP reporter gene expression. (B) Lentiviruses were produced in adherent HEK-293 cells grown in serum-free culture medium, using 15 µg DNA and 30 µl PEIpro® per 75 cm² flask. Virus yields were determined by titration of the supernatant 48 hr after transfection.

## 3. Efficient Virus Production in Any System at Any Scale

Figure 3: PEIpro® Is the Reagent of Choice for Virus Production Runs in Most Cell Culture Systems in Both Adherent and Suspension Cells, From Small Scale to Large Scale

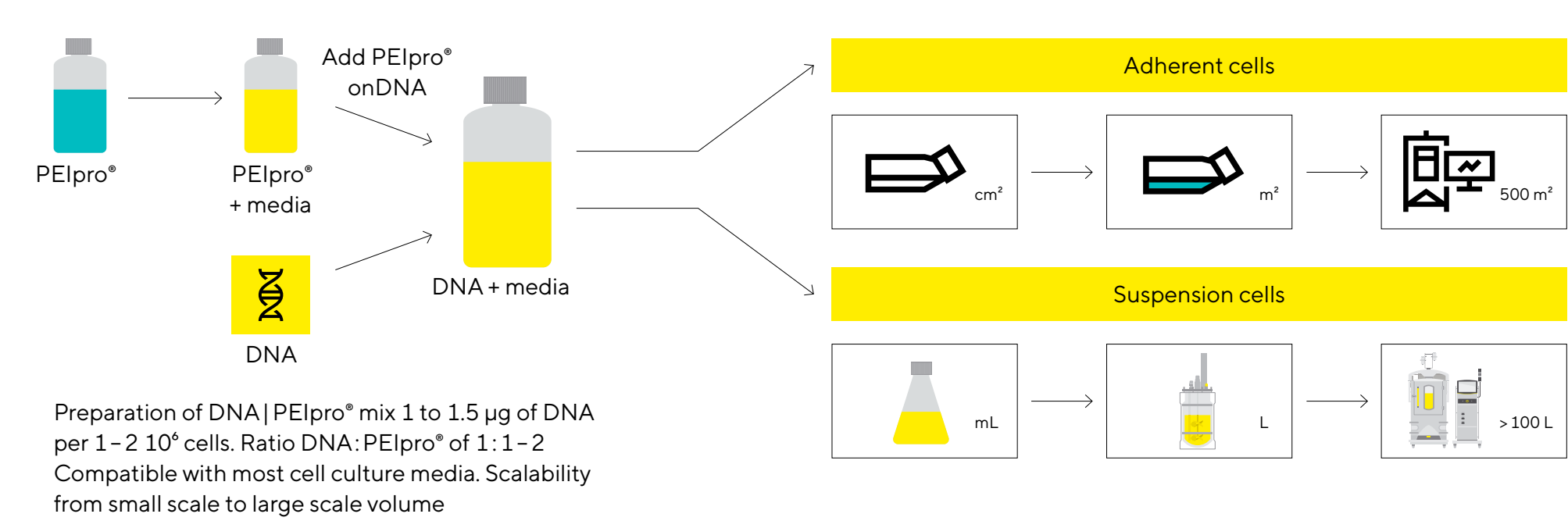
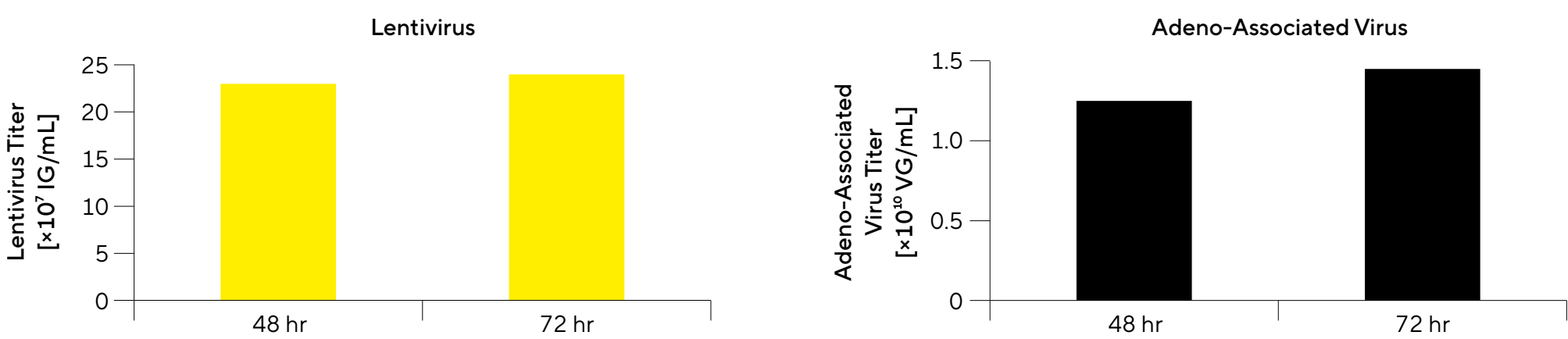


Figure 4: Lentivirus and AAV Production in HEK-293T and HEK-293 Cells Grown in Suspension in BalanCD® HEK-293 (Irvine Scientific®)



Note. HEK-293T (lentivirus) and HEK-293 (AAV) cells were thawed directly into each medium and passaged every 3 to 4 days before going into a 2 L benchtop bioreactor. Cells were seeded and cultured for 3 days before being transfected with PEIpro® (Polyplustransfection®). For transfection, four plasmids were used for lentivirus and three plasmids were used for AAV. Lentiviral and AAV titer were measured 48 and 72 hours post-transfection (Data kindly provided by Généthon).

Figure 5: PEIpro® to Simplify Scale-Up and to Ensure Reproducible Virus Production Yields in iCELLis® Nano Bioreactor

