SARTURIUS

PEIpro® Transfection Reagent

For Small- to Large-Scale Viral Vector Production



Product Information

PElpro® is the leading PEl-based DNA transfection reagent that offers flexibility and scalability to develop a robust and sustainable viral vector manufacturing process. PElpro® benefits from extensive research in PEl polymer chemistry and formulation to achieve the highest transfection efficiency in adherent and suspension cell culture systems. PElpro® allows a smooth transition from initial process development to large-scale manufacturing of therapeutic viral vectors with three quality grades available: PElpro®, PElpro®-HQ, and PElpro®-GMP. PElpro®-GMP is the highest quality grade PEl-based reagent on the market and is provided with full regulatory documentation to support drug approvals.

Features and Benefits

- Gold Standard: Optimized PEI-based transfection reagent for scalable viral vector production
- Reproducible: Equivalent viral vector production guaranteed through three quality grades
- **GMP-Grade:** Manufactured in compliance with ICH Q7 guidelines
- Residual Test: Proprietary residual test to mitigate risks and ensure patient safety

Introduction

Relevant Applications

- Viral vector production (AAV | LVV | others)
- Cell and gene therapy

Relevant Process Steps

Transfection (upstream process)

Three Quality Grades Support Seamless Transition From Process Development to Commercialization

The extensive research and development invested in PElpro® and the Polyplus (now part of Sartorius) expertise in the manufacturing process has made PElpro® the gold standard among commercially available PEls for cell and gene therapy. Three quality grades are available to guarantee the level of process qualification, quality controls, traceability, and documentation required for each project stage, enabling a seamless transition from process development to commercialization (Table 1). Both PElpro® and PElpro®-HQ are produced in an ISO 9001 environment, ensuring robust quality control. The PElpro®-GMP is produced in GMP-certified facilities following ICH Q7 guidelines that ensure a validated aseptic manufacturing process.

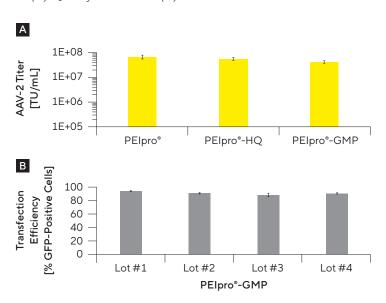
Table 1: Comparison of Three Quality Grades of PElpro® Transfection Reagent

Characteristics	PEIpro [®]	PElpro°-HQ	PElpro°-GMP	
Composition	Gold standard PEI-based transfection reagent for viral vector manufacturing used in cell and gene therapy. Ready-to-use, chemically-defined, and free of animal-derived components.			
Recommended use	Discovery to process development	Pre-clinical phase	Clinical trials and commercialization	
Quality grade	Research use only	Highly qualified grade	cGMP grade	
Packaging and size	 1.5 mL vial 10 mL bottle 100 mL bottle 1 L bottle 	100 mL bottle1 L bottle	 10 × 10 mL bottles 100 mL bottle 1 L bag (closed system) 	
Fill and finish manufacturing process	Filtration 0.22 μm	Filtration 0.22 μm	Sterile filtrationValidated aseptic processSimulations	
Quality controls	Standard QCs	Extended QCs to assess identity, potency, purity and safety	Validated QCs according to European pharmacopeia assessing identity, potency, purity and safety	
Product documentation	Certificate of analysisCertificate of originNon-hazardous product statement	Certificate of analysisCertificate of originNon-hazardous product statement	 Certificate of analysis Certificate of compliance Certificate of origin Non-hazardous product statement 	
Regulatory documentation	n/a	Batch production documentationQuality agreement	 Drug master file (DMF) Chemistry, manufacturing and control (CMC) section Protocols for identity, concentration and activity testing Quality agreement 	
Audit	According to ISO 9001 2015	According to ISO 9001 2015	According to ICH Q7 Eudralex Volume 4 Part II and Annex I	

A Scalable and Reproducible Transfection Reagent For a Robust Manufacturing Process

To ensure robust manufacturing, scaled-up or scaled-out therapeutic virus production with PElpro®-HQ or PElpro®-GMP is designed to maintain the production yields achieved with R&D grade PElpro® during process development. PElpro® is manufactured and formulated using a highly controlled and strict manufacturing process. By ensuring reproducibility from one run to the next, PElpro® greatly facilitates process standardization, a key feature of reliable large-scale production. Moreover, PElpro® is released using advanced quality controls, including a specification for transfection efficiency that supports excellent lot-to-lot consistency (Figure 1).

Figure 1: High Reproducibility with PElpro® Regardless of (A) Quality Grade and (B) Production Lot

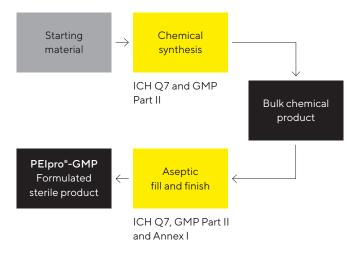


Note. (A) Suspension HEK-293T cells were seeded at 1×10^6 cells/mL in a commercial culture medium and transfected with either PElpro®, PElpro®-HQ or PElpro®-GMP reagents following the same protocol for each product. AAV-2 titers were measured 72 hours after transfection using a GFP reporter gene expression. (B) Suspension HEK-293T cells were seeded at 1×10^6 cells/mL in a commercial culture medium and transfected with four different lots of PElpro®-GMP with a GFP-expressing plasmid. Transfection efficiency was measured 48 hours post-transfection by flow cytometry.

Risk Mitigation With PElpro®-GMP

Introducing any starting or raw | ancillary material in the manufacturing process of cell and gene therapies adds potential risk factors. Using PEIpro®-GMP as the transfection reagent to produce viral vectors increases confidence in the quality and safety of the process and product. PEIpro®-GMP is manufactured in compliance with international GMP guidelines ICH Q7 and correspondingly with EU "Guidelines for good manufacturing practices for medicinal products for human and veterinary use" (ICH Q7 and EudraLex Vol 4, Part II and Annex I: Manufacture of Sterile Medicinal Products). PEIpro®-GMP 1 L is provided in Flexsafe® bags with MPC connectors and weldable tubing suitable for closed systems to decrease contamination risks. A residual test to analyze the presence of the transfection reagent in the drug product is available through selected service providers.

Figure 2: PElpro®-GMP manufacturing process



Note. PElpro®-GMP is provided with a full set of tools to ensure traceability with dedicated regulatory documentation and analytics with different protocols to support both IND and BLA submissions.

Technical Specifications

PEIpro®

Attribute	PElpro [®] 1.5 mL	PElpro [®] 10 mL	PElpro® 100 mL	PElpro® 1 L
Quality grade	Research grade	Research grade	Research grade	Research grade
Type of container	Vial	Bottle	Bottle	Bottle
Material	Polypropylene	Polypropylene	Polypropylene	Polypropylene
Neck diameter [mm]	n/a	13.7 ± 0.05	17.5 ± 0.05	27.4 ± 0.08
Height with closure [mm]	n/a	58.4 ± 1.5	100.8 ± 2.0	215.9 ± 2.5
Bottle diameter [mm]	n/a	24.9 ± 1.0	50.3 ± 1.5	91.4 ± 2.0

PEIpro®-HQ

Attribute	PEIpro°-HQ 100 mL	PElpro°-HQ 1 L
Quality grade	Highly qualified	Highly qualified
Type of container	Bottle	Bottle
Material	Polypropylene	Polypropylene
Neck diameter [mm]	17.5 ± 0.05	27.4 ± 0.08
Height with closure [mm]	100.8 ± 2.0	215.9 ± 2.5
Bottle diameter [mm]	50.3 ± 1.5	91.4 ± 2.0

PEIpro®-GMP

Attribute	PEIpro°-GMP 10 × 10 mL	PEIpro°-GMP 100 mL	PEIpro®-GMP 1 L	PEIpro®-GMP Satellite Bag 100 mL
Quality grade	GMP	GMP	GMP	GMP (for ID test)
Type of container	Bottle (×10)	Bottle	Flexsafe® 2D Bag 1 L	Flexsafe® 2D Bag 1 L
Material	PETG (sterile)	PETG (sterile)	S80 Polyethylene film	S80 Polyethylene film
Neck diameter [mm]	14	28	n/a	n/a
Height [mm]	64	110	n/a	n/a
Bottle diameter [mm]	38	54	n/a	n/a
Bag dimension (W × H) [mm]	n/a	n/a	240 × 310	240 × 310
Connector	n/a	n/a	MPC (coupling male 1:4")	MPC (coupling male 1:4")
Other	n/a	n/a	Weldable tubing	Weldable tubing

Ordering Information

Item	Volume	Package	Order Number
PElpro®	1.5 mL	Vial	101000017
PElpro®	10 mL	Bottle	101000033
PElpro®	100 mL	Bottle	101000026
PElpro®	1 L	Bottle	101000029
PEIpro®-HQ	100 mL	Bottle	101000052
PEIpro®-HQ	1 L	Bottle	101000039
PEIpro®-GMP	10 × 10 mL	Bottle	102000131
PEIpro®-GMP	100 mL	Bottle	102000121
PEIpro®-GMP	1 L	Bag	102000008
PEIpro®-GMP satellite bag	100 mL	Bag	102000004

Germany

Sartorius Stedim Biotech GmbH August-Spindler-Strasse 11 37079 Goettingen Phone +49 551 308 0

USA

Sartorius Stedim North America Inc. 565 Johnson Avenue Bohemia, NY 11716 Toll-Free +1 800 368 7178

France

Polyplus®—Now part of Sartorius 75 Rue Marguerite Perey 67400 Illkirch Phone +33 390 406 180

For more information, visit

sartorius.com/transfection-reagents-plasmids polyplus-sartorius.com/products/peipro