

FectoVIR®-LV Transfection Reagent

Lentiviral Vector
Manufacturing



Product Information

FectoVIR®-LV is an animal component-free, next-generation transfection reagent designed to enhance the production of lentivirus (LV) vectors in HEK-293 cell systems. Ideal for large-scale manufacturing, FectoVIR®-LV enables a reduction of the complexation volume and increases complex stability. These benefits allow manufacturers to increase the number of doses produced per batch while decreasing LV manufacturing costs, increasing patient accessibility. FectoVIR®-LV GMP and a residual test are currently under development to support GMP applications.

Features and Benefits

- Productivity: Maximize LV vector titers
- Cost-effectiveness: Reduce your cost per batch
- Scalability: Manufacture at large scale with low complexation volume and long complex stability
- Time-saving: Use DOE to help you optimize your process in a record time

Introduction

Relevant Applications

- Cell and gene therapy
- Gene-modified cell therapy
- CAR-T, CAR-NK, and CAR-TCR
- Gene-modified stem cell therapy
- LV vector production
- Academic research using LV vectors

Relevant Process Steps

- Upstream process
- Transfection

Figure 1: *FectoVIR®-LV Demonstrates High Productivity for LV Vector Manufacturing Compared to Competitor Reagents*

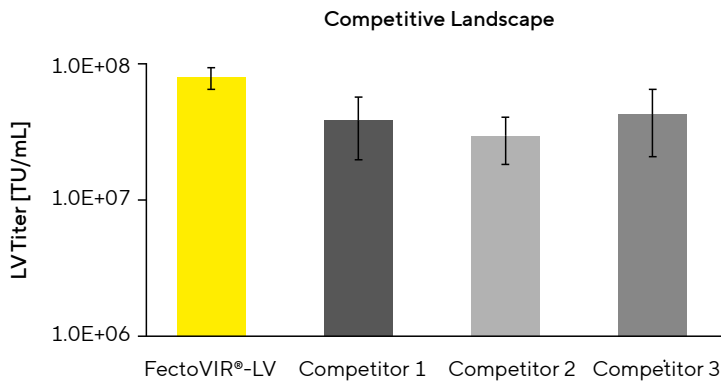


Figure 2: *FectoVIR®-LV Enables the Use of Lower DNA Quantities While Optimizing Titers to Improve Process Economics*

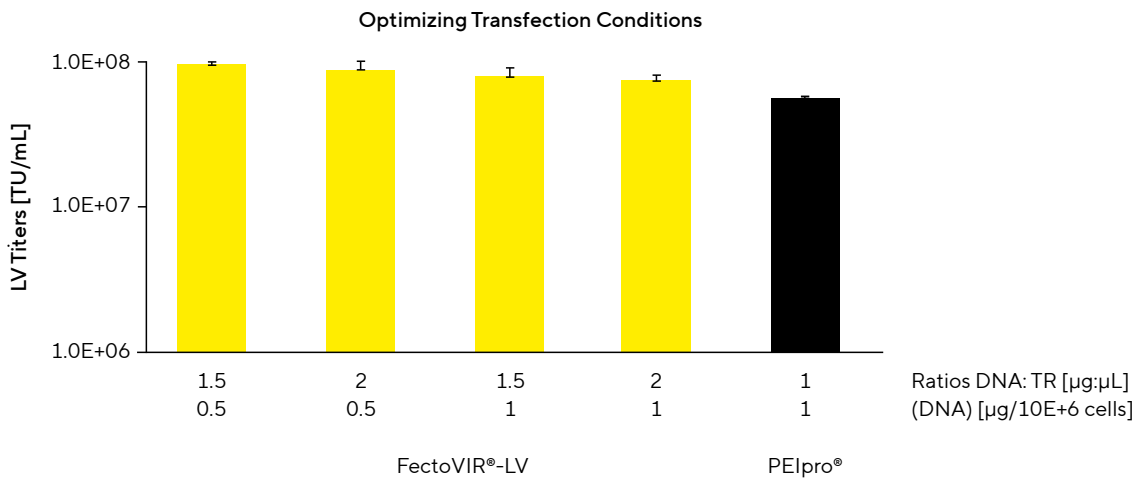


Figure 3: FectoVIR®-LV Offers Robust Scalability by Reducing the Complexation Volume and Increasing the Complex Stability

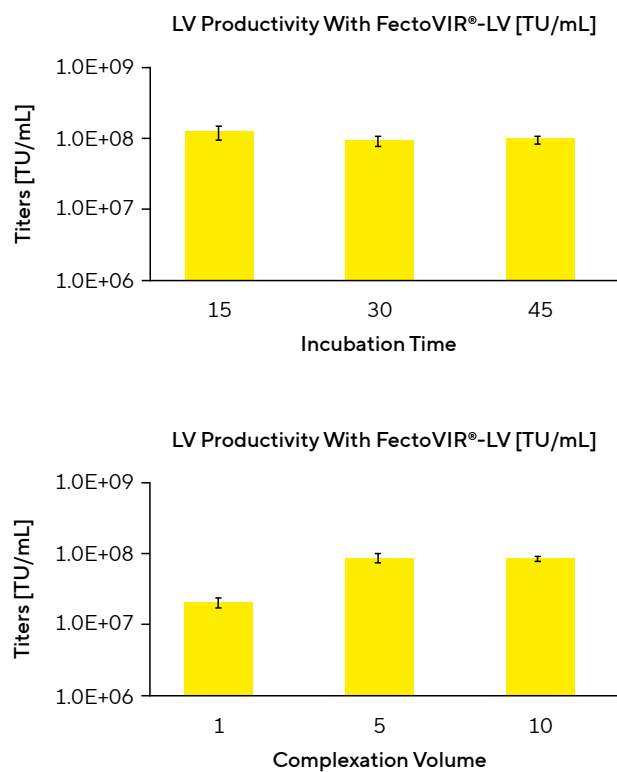
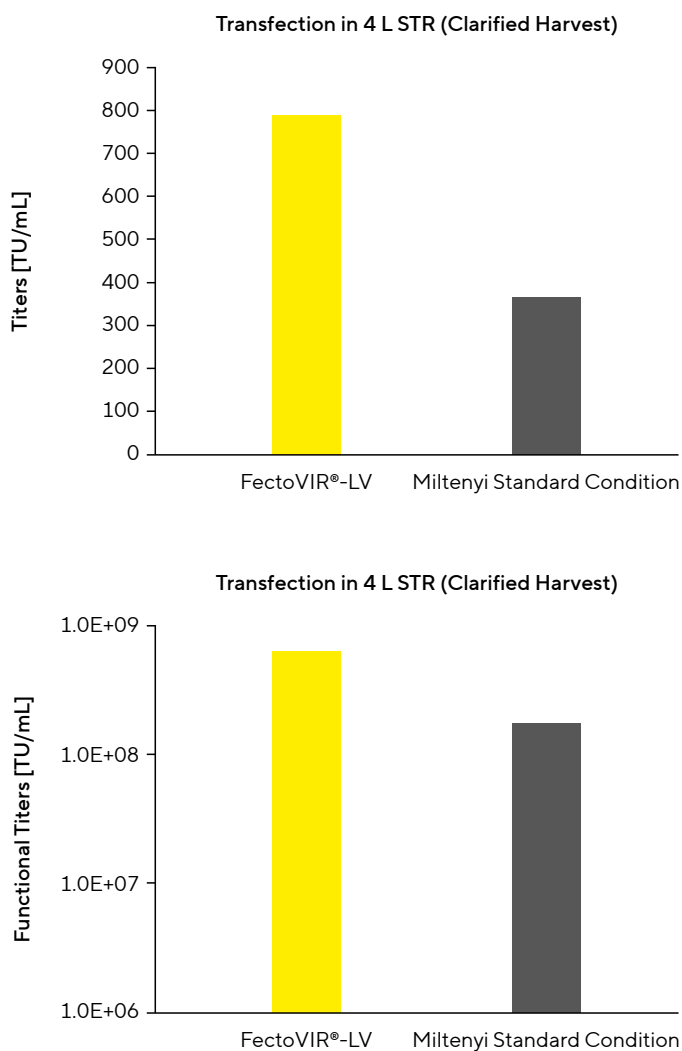


Figure 4: FectoVIR®-LV Confirmation Run in a 4 L Flask After DOE Optimization



Technical Specifications

Attribute	FectoVIR®-LV 1 mL	FectoVIR®-LV 10 mL	FectoVIR®-LV 100 mL
Quality grade	Research grade	Research grade	Research grade
Transfected cell culture volume*	500 mL	5 L	50 L
Type of container	Vial	Bottle	Bottle
Material	Polypropylene	Polypropylene	Polypropylene
Neck diameter [mm]	n/a	13.7 ± 0.05	17.5 ± 0.05
Height with closure [mm]	n/a	58.4 ± 1.5	100.8 ± 2.0
Height without closure [mm]	n/a	55.9 ± 1.5	98.6 ± 2.0
Bottle diameter [mm]	n/a	24.9 ± 1.0	50.3 ± 1.5

* Average volume of cell culture to be transfected may vary depending on the transfection conditions.

Ordering Information

Item	Description	Volume	Order Number
FectoVIR®-LV 1 mL	Vial	1 mL	101000187
FectoVIR®-LV 10 mL	Bottle	10 mL	101000188
FectoVIR®-LV 100 mL	Bottle	100 mL	101000189

 **To Order Polyplus® Products, Visit**
polyplus-sartorius.com
polyplus-sartorius.com/products/fectovir-lv

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 3 90 40 61 80

 **For more information, visit**
sartorius.com
polyplus-sartorius.com
polyplus-sartorius.com/products/fectovir-lv

Specifications subject to change without notice.

©2024 Sartorius Stedim Biotech GmbH, August-Spindler-Strasse 11, 37079 Goettingen, Germany

Status: 03 | 2024