

# pPLUS® AAV-RC Range

Ready-to-use RepCap  
Plasmid for AAV Production



## Product Information

pPLUS® AAV-RC are ready-to-use RepCap plasmids designed for production of various AAV serotypes through transient transfection of HEK293 cells in suspension. It has been optimized for use with pPLUS® AAV-Helper and FectoVIR®-AAV and can be used with various HEK293 cell lines.

## Features and Benefits

- **Enhanced AAV Productivity When Used With pPLUS® AAV-Helper:** Enhancement of VG titers and TU titers
- **Readily Available:** Ready-to-use product, available off-the-shelf in three pack sizes from 1 to 100 mg
- **Royalty-Free Even for Commercial Use**

# Introduction

## Relevant Applications

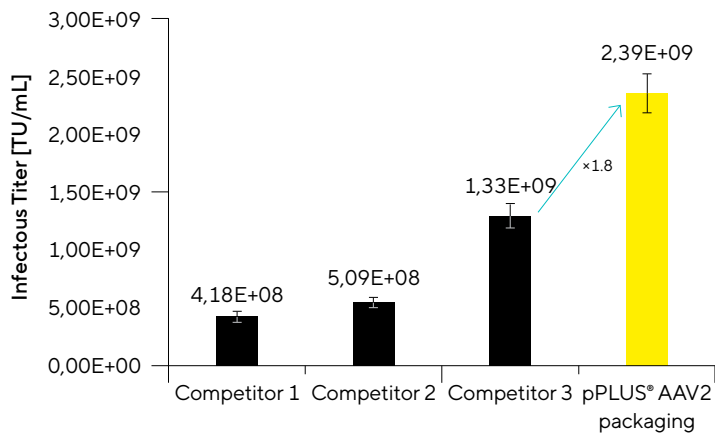
- Cell and Gene Therapy
- AAV production
- Academic Research

## Relevant Process Steps

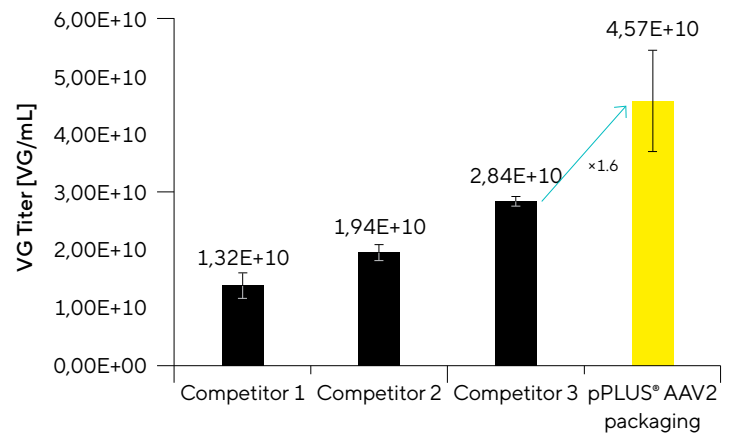
- Upstream Process
- Transfection

**Figure 1:** The Use of pPLUS® AAV Packaging System for AAV2 leads to 1.8 to 5.7 fold increase in TU and 1.6 to 3.5 fold increase in VG

Performance of pPLUS® AAV Packaging Plasmids AAV2 Infectivity

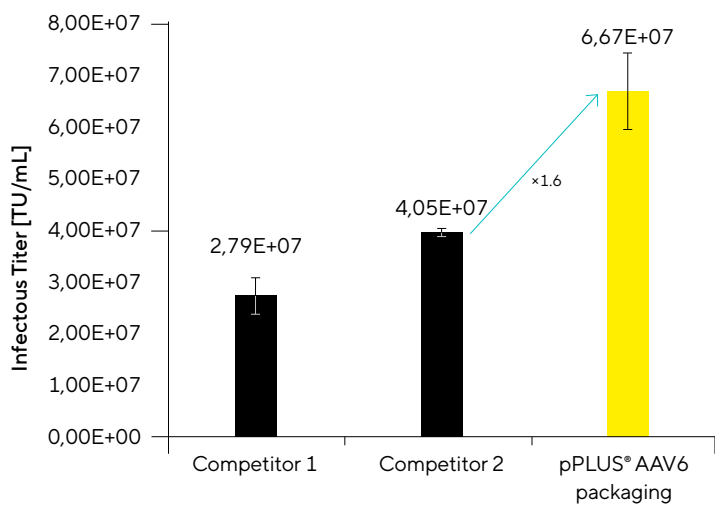


Performance of pPLUS® AAV Packaging Plasmids AAV2 Genomic Titer

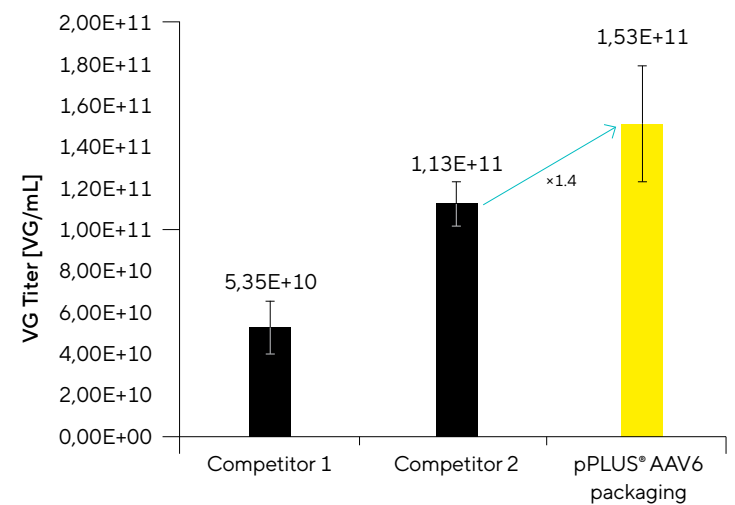


**Figure 2:** The Use of pPLUS® AAV Packaging System for AAV6 leads to 1.6 to 2.4 fold increase in TU and 1.4 to 2.9 fold increase in VG

Performance of pPLUS® AAV Packaging Plasmids AAV6 infectivity



Performance of pPLUS® AAV Packaging Plasmids AAV6 Genomic Titer



Note. AAV were produced using pPLUS® AAV packaging plasmids (Helper and RepCap) in optimized condition following our DOE experiment vs competitors packaging plasmids and their recommended condition. The plasmids GFP transgene used were the same for all experiments and were internally engineered. AAV were internally produced in VPC2.0 cells, VPM medium (ThermoFisher) with FectoVIR®-AAV. VG titers were measured through dPCR and infectious titers were measured after transduction of HT-1080 cells.

# Technical Specifications

Serotype	AAV2, AAV5, AAV6
Plasmid size	6100 – 6300 bp
Antibiotic resistance gene	Kan/neoR
Appearance	Liquid
Storage temperature	-20°C

## Ordering Information

Item	Description	Volume	Order Number
pPLUS® AAV-RC2 1 mg	RepCap2 plasmid for AAV production	1 mg vial	101000214
pPLUS® AAV-RC2 10 mg	RepCap2 plasmid for AAV production	10 mg vial	101000215
pPLUS® AAV-RC2 100 mg	RepCap2 plasmid for AAV production	100 mg vial	101000216
pPLUS® AAV-RC5 1 mg	RepCap5 plasmid for AAV production	1 mg vial	101000242
pPLUS® AAV-RC5 10 mg	RepCap5 plasmid for AAV production	10 mg vial	101000243
pPLUS® AAV-RC5 100 mg	RepCap5 plasmid for AAV production	100 mg vial	101000244
pPLUS® AAV-RC6 1 mg	RepCap6 plasmid for AAV production	1 mg vial	101000245
pPLUS® AAV-RC6 10 mg	RepCap6 plasmid for AAV production	10 mg vial	101000246
pPLUS® AAV-RC6 100 mg	RepCap6 plasmid for AAV production	100 mg vial	101000247

**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](https://sartorius.com/transfection-reagents-plasmids)