

# pPLUS<sup>®</sup> AAV-RC2

Ready-to-use RepCap  
Plasmid for AAV2 Production



## Product Information

pPLUS<sup>®</sup> AAV-RC2 is a ready-to-use RepCap plasmid designed for production of AAV2 through transient transfection of HEK293 cells in suspension. It has been optimized for use with pPLUS<sup>®</sup> AAV-Helper and FectoVIR<sup>®</sup>-AAV and can be used with various HEK293 cell lines.

## Features and Benefits

- **Enhanced AAV Productivity When Used With pPLUS<sup>®</sup> 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

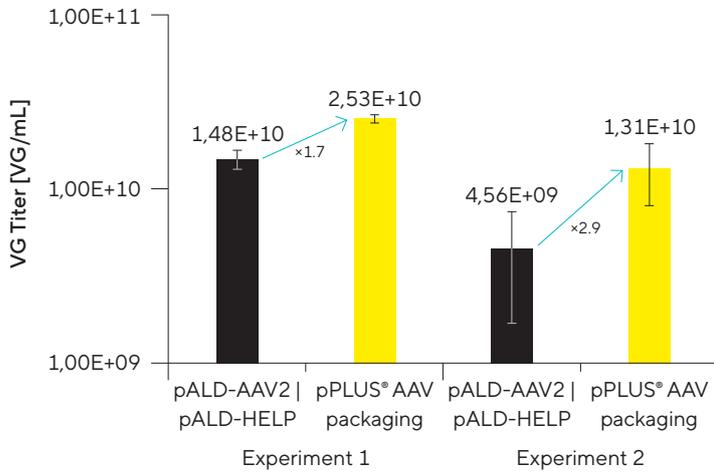
- Cell and Gene Therapy
- AAV production
- Academic Research

## Relevant Process Steps

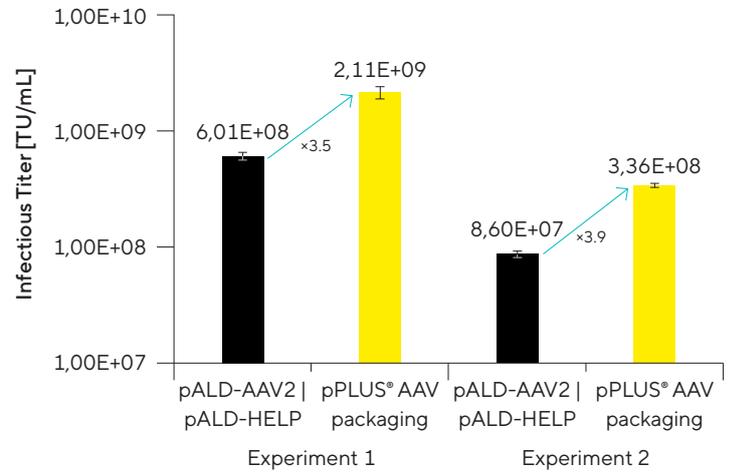
- Upstream Process
- Transfection

**Figure 1:** The Use of pPLUS® AAV Packaging System Leads to A 1.7-to-2.9-Fold Increase in VG Tires and 3.5 to 3.9 Fold Increase in TU Titors vs the Competitor Control

Performance of pPLUS® AAV Packaging Plasmids AAV2 Genomic Titer



Performance of pPLUS® AAV Packaging Plasmids AAV2 Infectivity



Note. AAV2 were produced using pPLUS® packaging plasmids vs pALD (Aldevron) packaging plasmids and an internally engineered pGFP in VPC2.0 cells | VPM medium (ThermoFisher) with FectoVIR®-AAV. VG titers were measured through qPCR and infectious titers were measured after transduction of HT-1080 cells.

# Technical Specifications

Serotype	AAV2
Plasmid size	6249 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

**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)