

# pPLUS® AAV-Helper, novel engineered pHelper plasmid to improve yield and quality of several AAV serotypes in suspension cell culture systems.

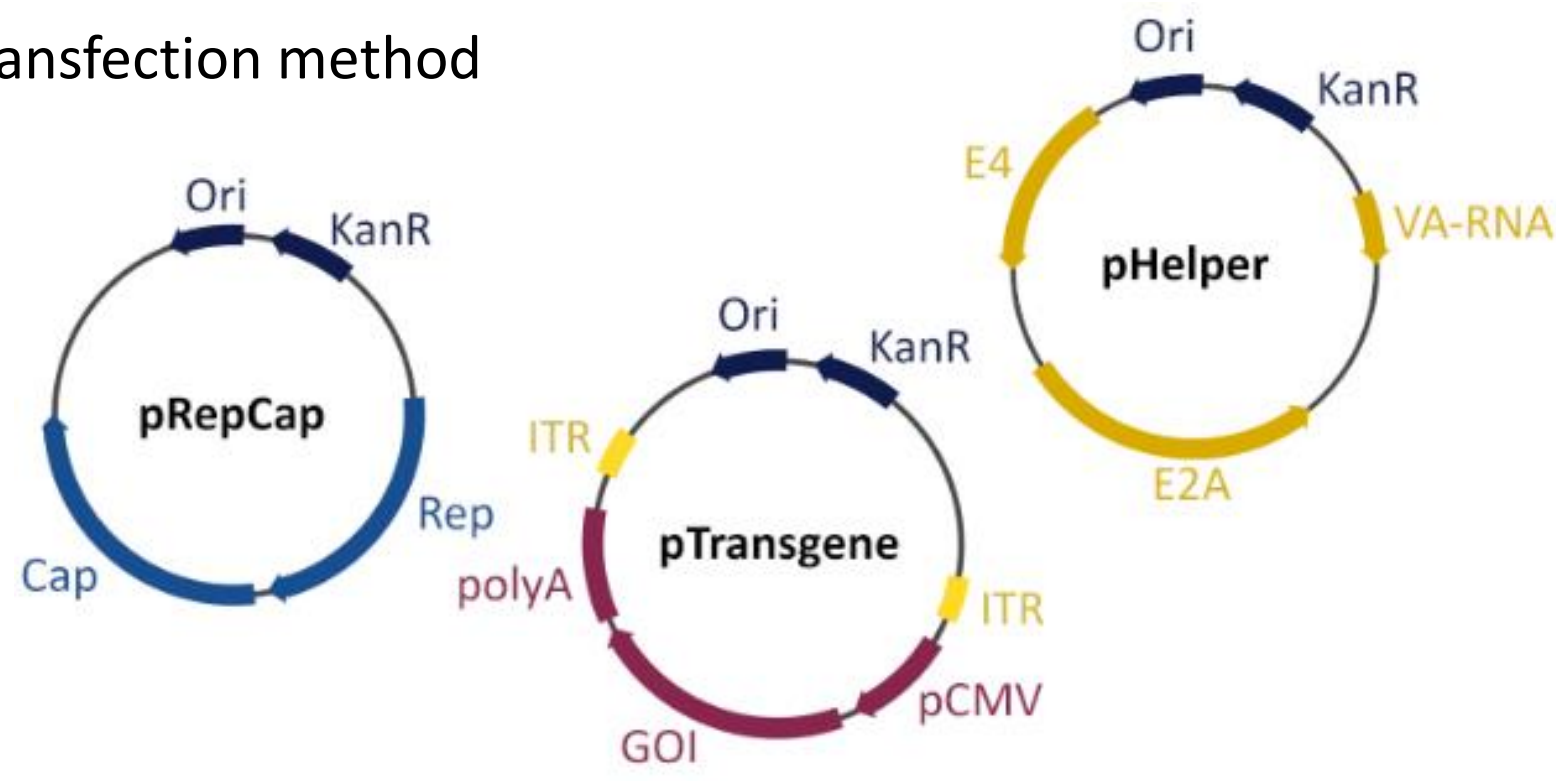
Eric Mauro<sup>2</sup>, Jonathan Havard<sup>1</sup>, Laure Robert<sup>1</sup>, Alengo Nyamay'antu<sup>1</sup>, Carine Morel<sup>2</sup>, Julie Regnery<sup>3</sup>, Elisa Berera<sup>3</sup>, Samantha Convers<sup>3</sup>, Pauline Schorter<sup>3</sup>, Quentin Bazot<sup>3</sup>, Sylvain Julien<sup>2</sup>, Patrick Erbacher<sup>1</sup>

<sup>1</sup> Polyplus®, Illkirch, France <sup>2</sup> Polyplus®, Loos, France <sup>3</sup> ABL, Lyon, France

Harnessing rAAVs as viral vectors for therapeutic transgene delivery still requires improvements in yields and specificity to lower vector doses, and therefore manufacturing cost, as well as to improve patient safety. To this end, our research is focused on developing novel technologies to ensure manufacturing of high yielding rAAV particles using transient transfection, as well as enhancing features of rAAV vectors that act on the overall size of packaged material and specificity of delivery. Here we present our state-of-the-art approach to design new helper plasmids (pHelpers) with the aim of improving both the infectivity (TU/mL) and the quality (full/empty ratio) of the viral particle obtained from suspension cultures. We took the opportunity to exploit our proprietary DNA assembly method technology to explore the synergies of multiple genetic features modularly assembled in synthetic plasmids. Comparison of the biological activity of several versions of rationally designed pHelpers led us to identify the optimal configuration able to outperform existing helper plasmids in every tested bioproduction conditions. Our expertise in DNA plasmid design and assembly together with our scalable transfection solutions for rAAV manufacturing gives us the potential to improve both productivity and specificity of gene therapy products.

## Introduction: rAAV production

### Triple transfection method



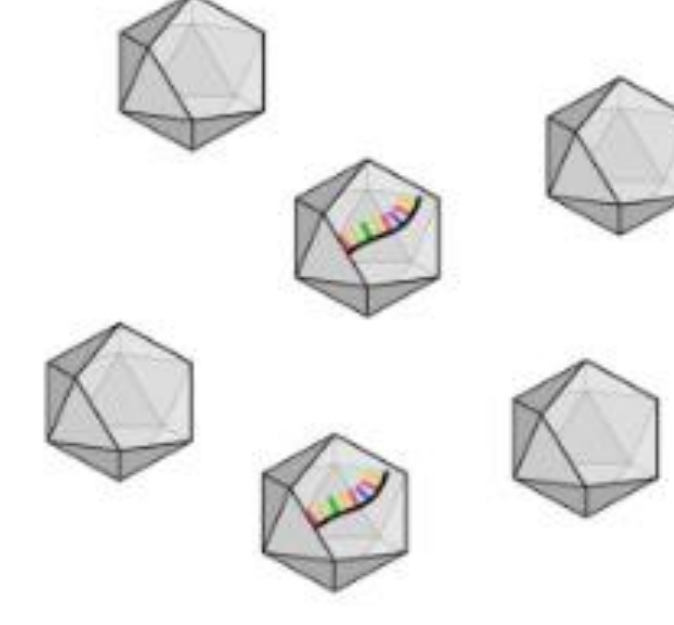
pHelper, pRepCap and pTransgene are transfected into HEK-293T cells

### Transfection



72h post-transfection, cells are lysed and viruses harvested

### Virus harvest



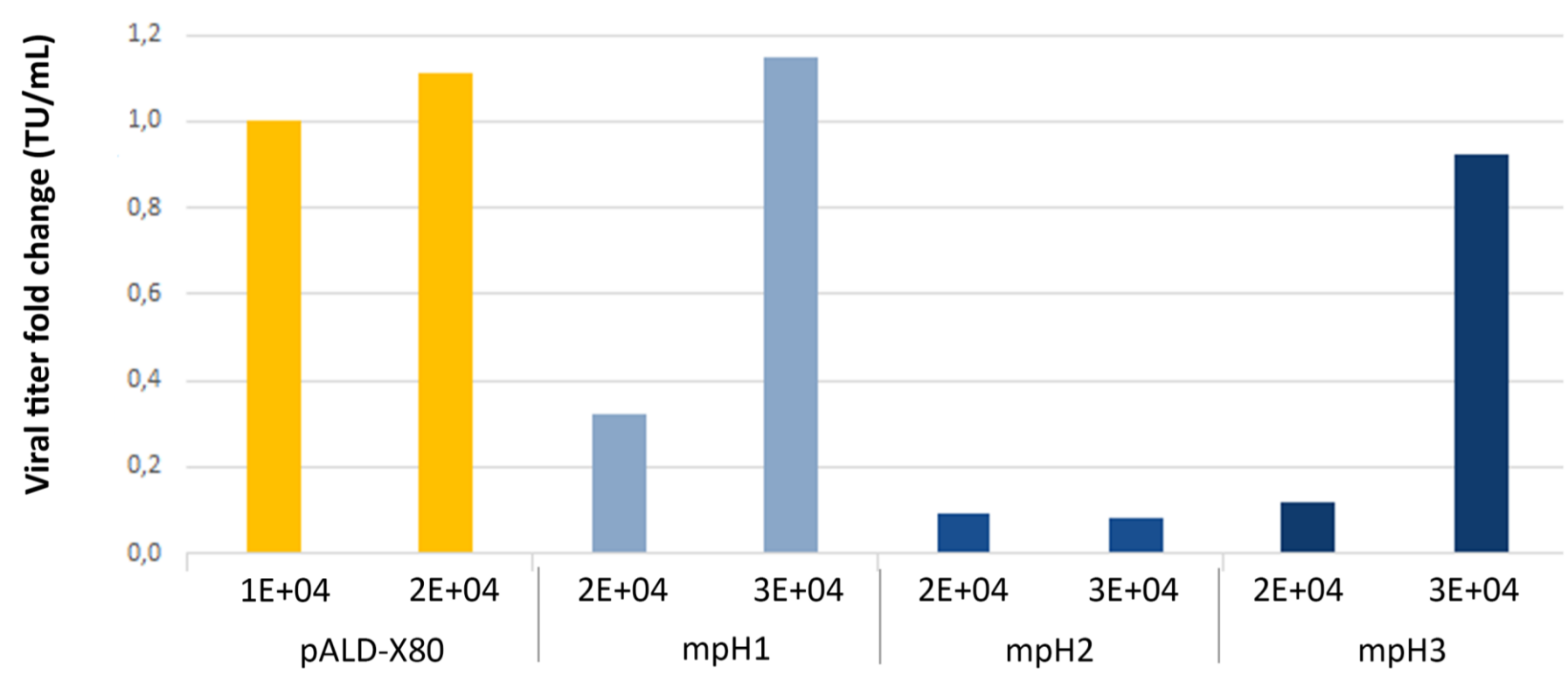
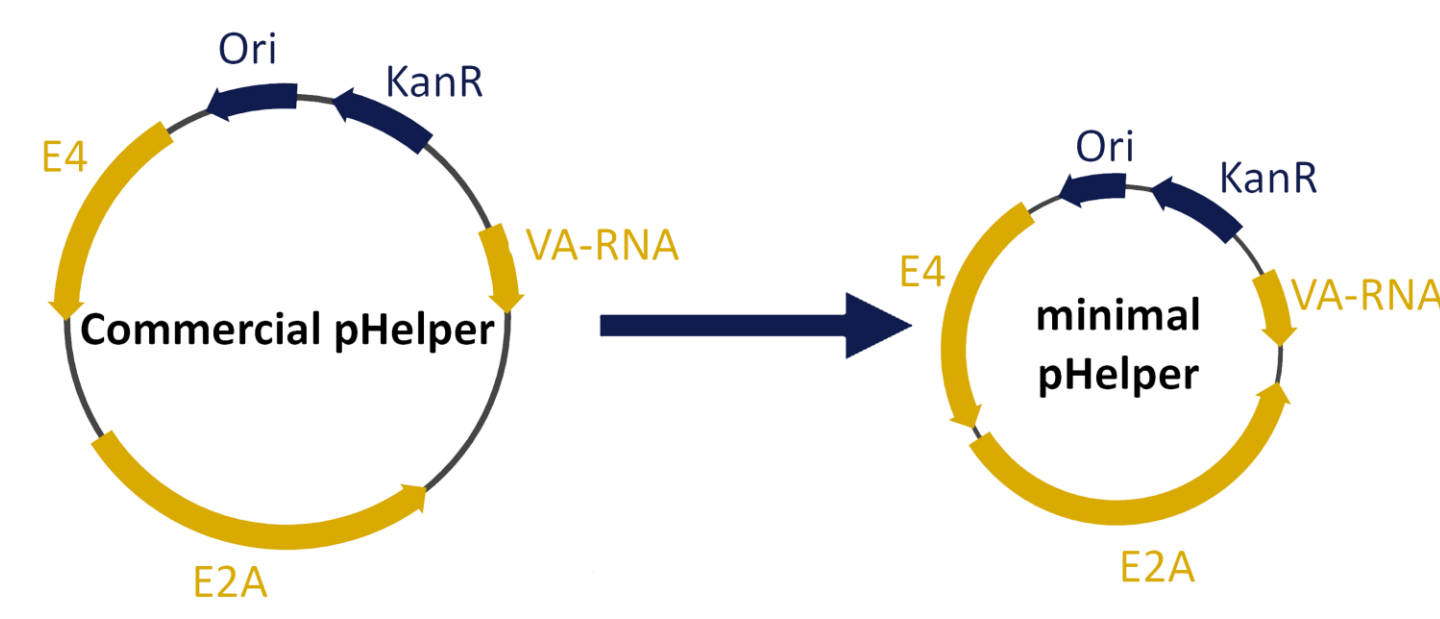
Viral titer is quantified by flow cytometry

### Strategy

- Various constructions of pHelper were made using e-Zyvec® plasmid assembly technology.
- Plasmids were transfected into HEK-293T cells and harvested 72h post-transfection.
- Viral titers (TU/mL) of the produced rAAV were then compared to identify the best pHelper for optimal rAAV production.

## Removing non essential elements

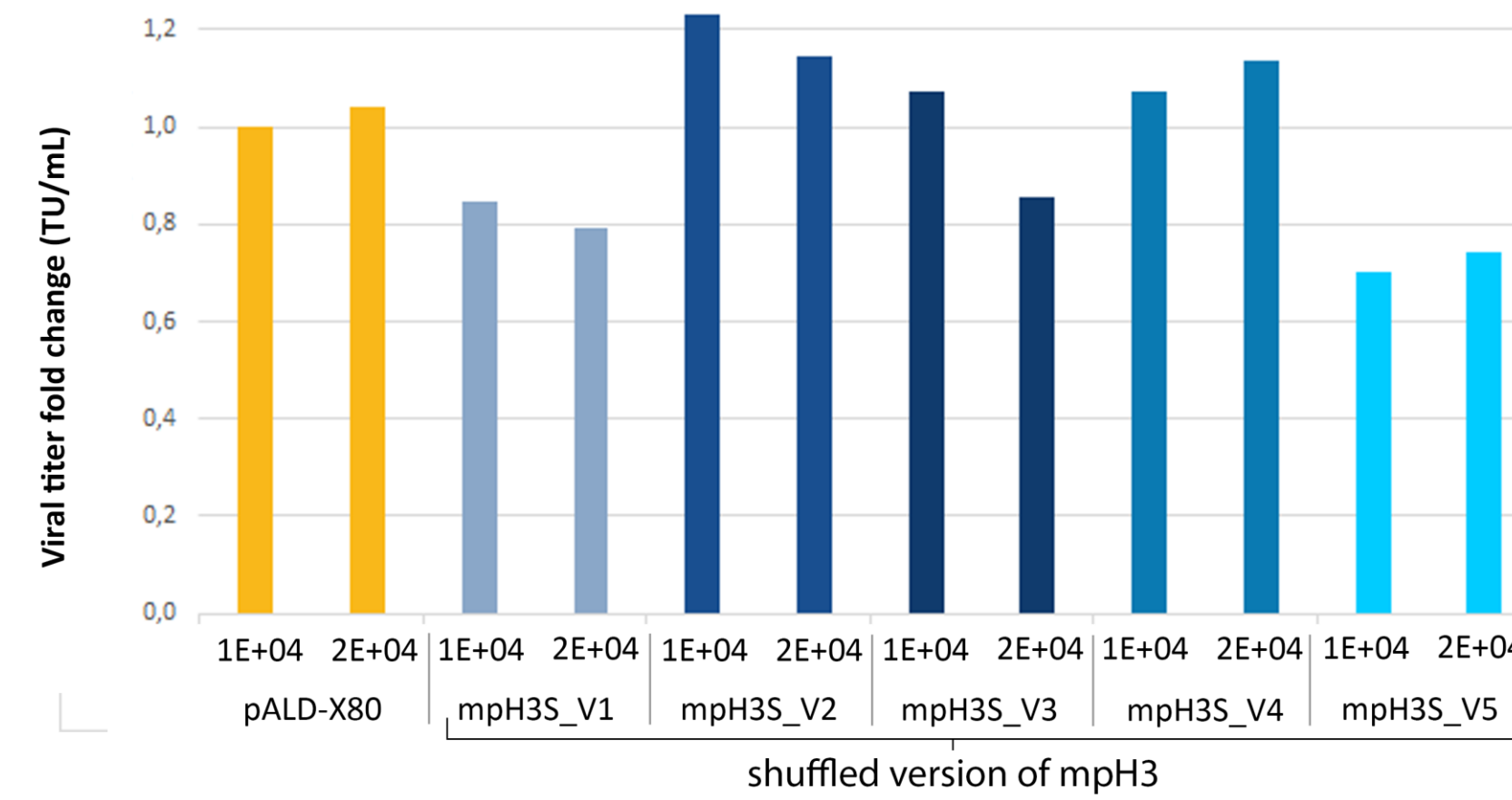
Construction of a minimal pHelper (mpH).



**Fig1:** Viral titers of rAAV2 (normalized to pALD-X80). Copy/cell of the pHelpers are shown above plasmids name.  
19 constructions tested, 3 series

## Shuffling

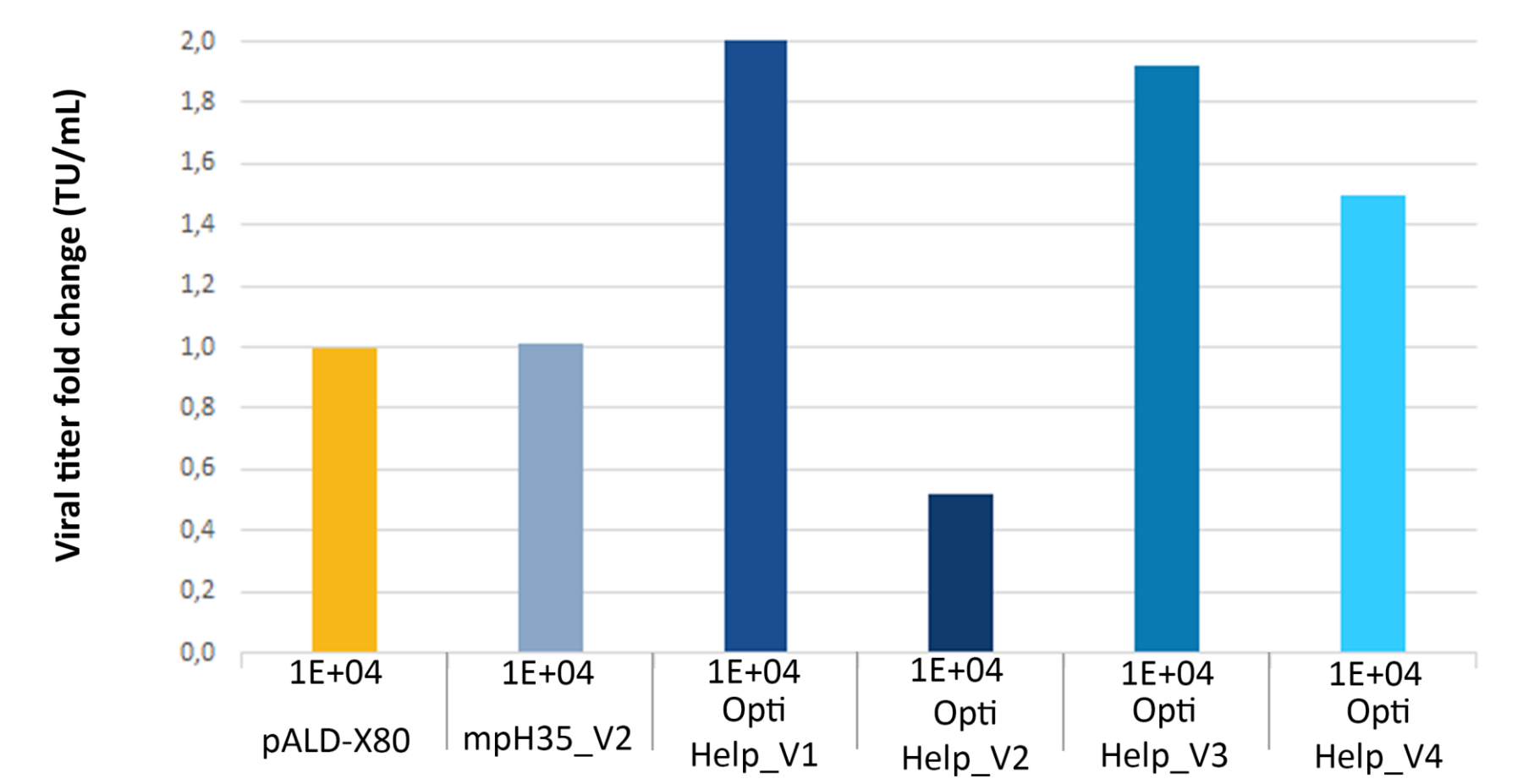
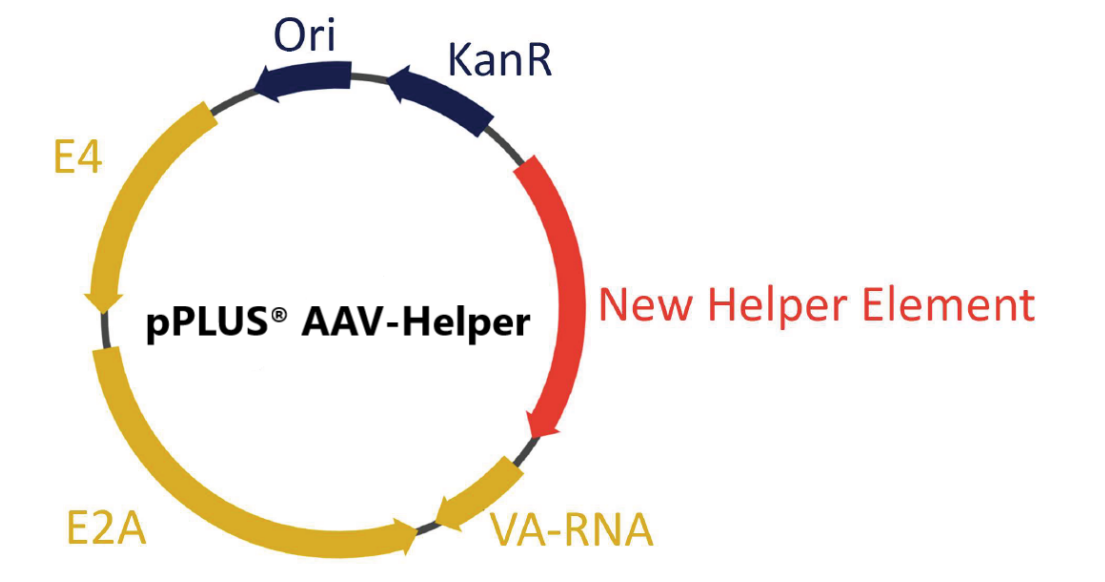
Re-organization of plasmid configuration to optimize its efficiency to produce rAAV.



**Fig2:** Viral titers of rAAV2 (normalized to pALD-X80). Copy/cell of the pHelpers are shown above plasmids name.  
9 constructions tested, 2 series

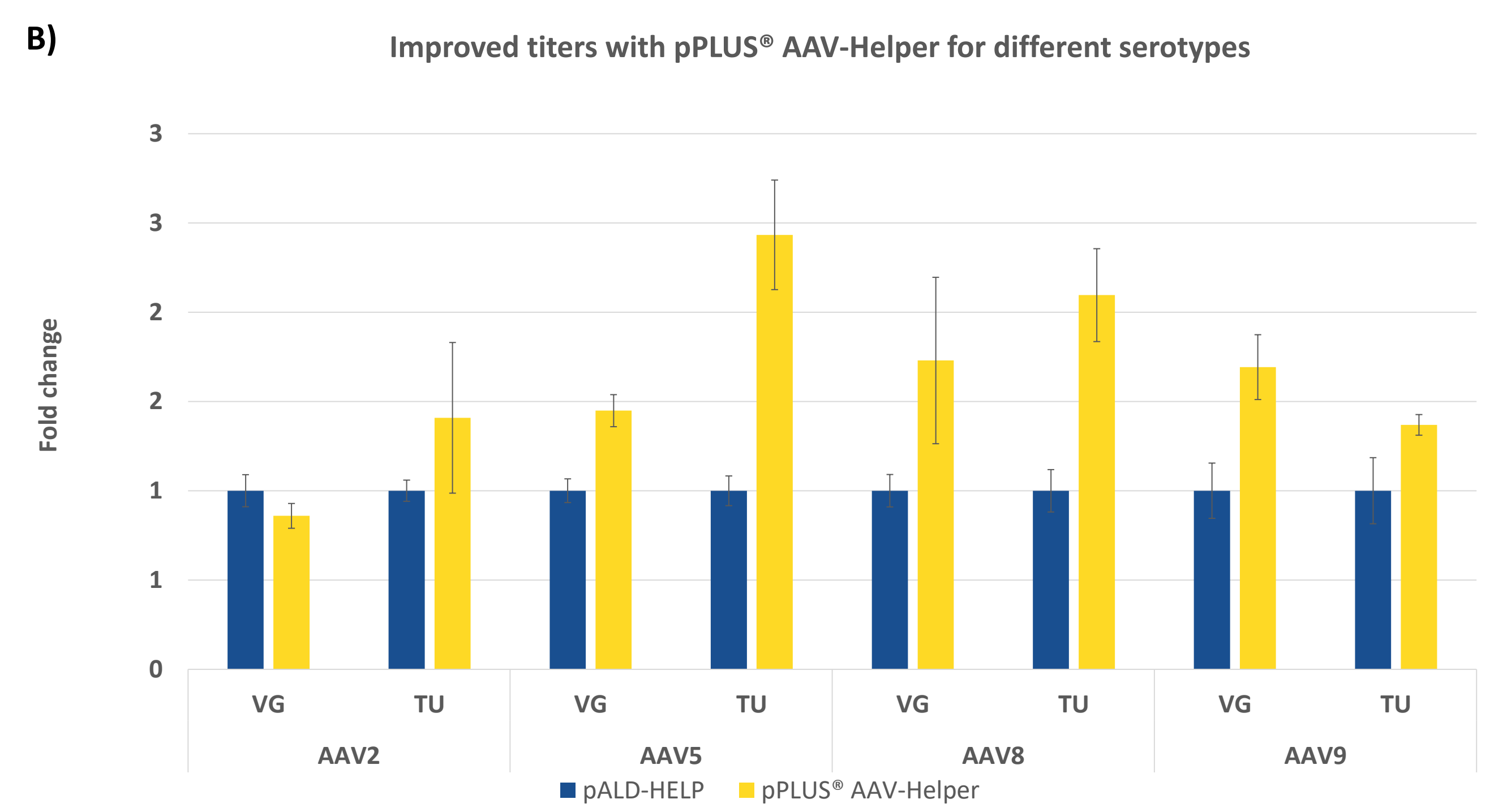
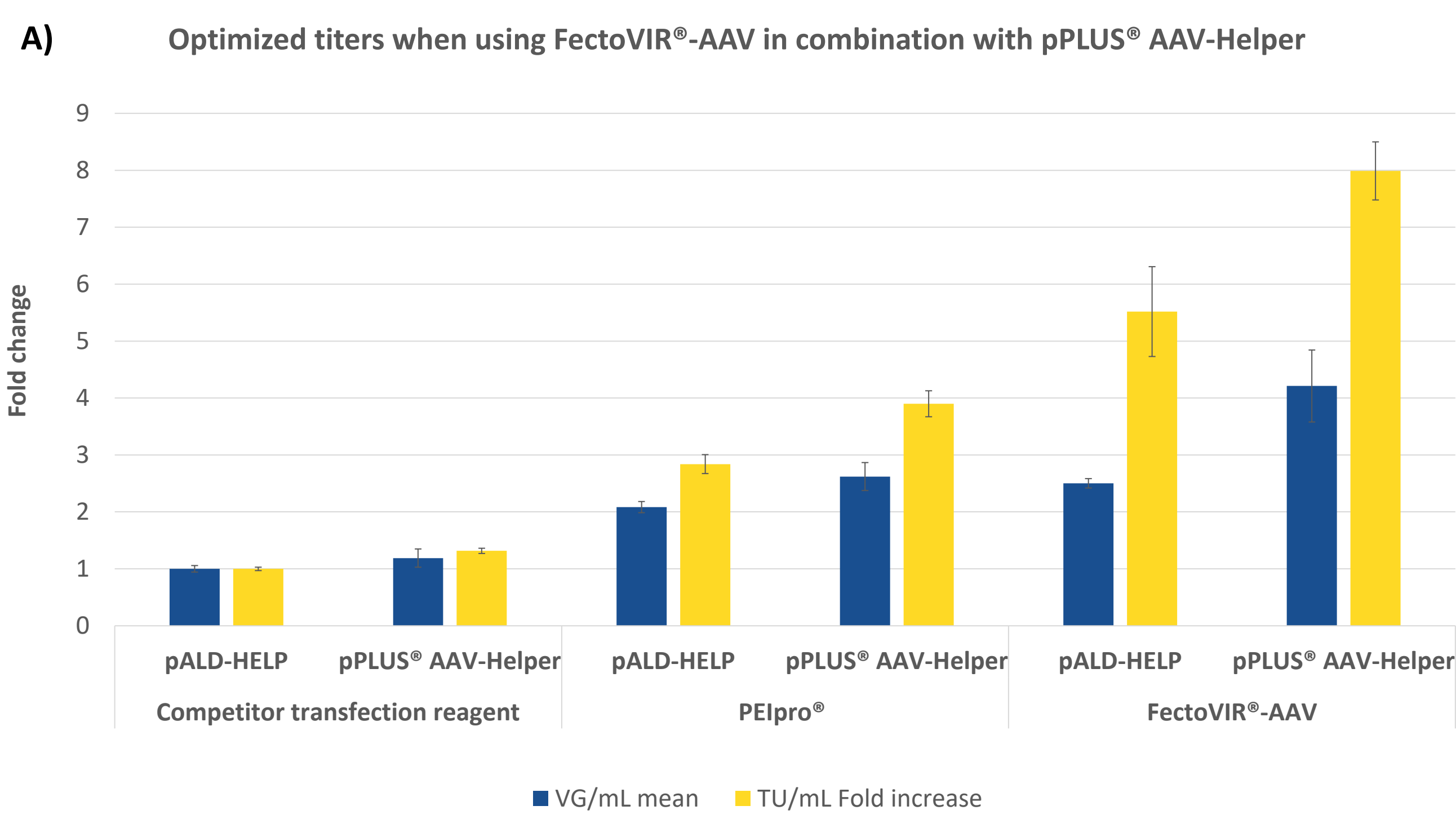
## Addition of new Helper elements

Identification of new sequences to add to the pHelper to boost rAAV production.

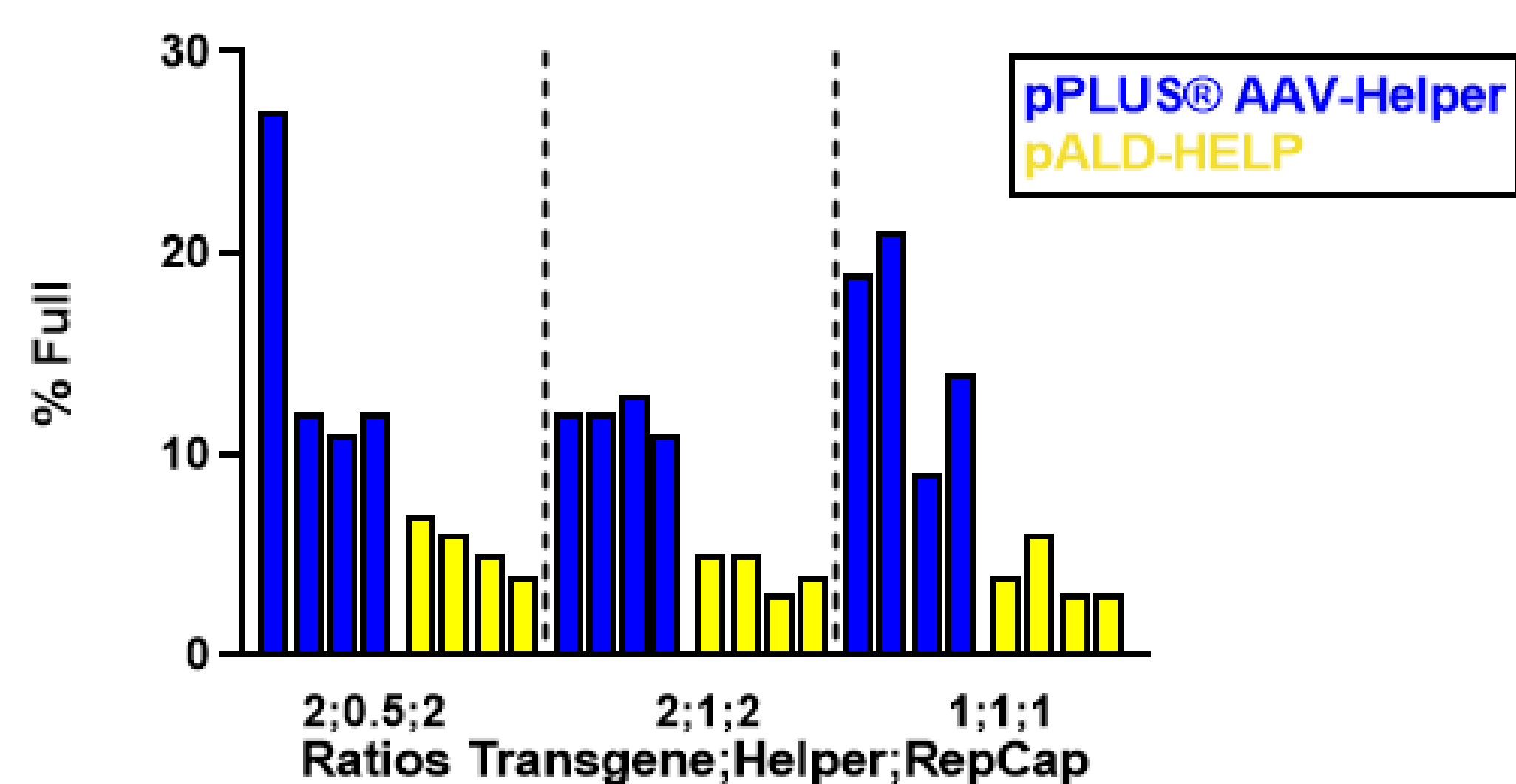


**Fig3:** Viral titers of rAAV2 (normalized to pALD-X80). Copy/cell of the pHelpers are shown above plasmids name.  
29 constructions tested, 5 series

## pPLUS® AAV-Helper to improve yield and quality of different AAV serotypes



**C) Improved full/empty particle ratio with pPLUS® AAV-Helper**



**Fig 4:** pPLUS® AAV-Helper is a novel helper plasmid for AAV production in suspension HEK-293 cell lines to improve production yields of all AAV serotypes tested (AAV2, AAV5, AAV8, AAV9).

**A) Improved productivity and infectivity with an optimized efficiency when used in synergy with FectoVIR®-AAV transfection reagent.** AAV9 were produced in HEK293T cells adapted in suspension in F17 medium and transfected with FectoVIR®-AAV. Transduction assay was performed on HEK293T cells.

**B) Improved titers with pPLUS® AAV-Helper for different serotypes.** AAV2, AAV5, AAV8 and AAV9 were produced in HEK293T cells adapted in suspension in F17 medium and transfected with FectoVIR®-AAV. Transduction assay was performed on HT-1080 cells. Improvement of titers are shown as fold increase in reference to pALD-HELP.

**C) Improved full/empty particle ratio with pPLUS® AAV-Helper.** AAV2 was produced in HEK293 suspension cells and transfected with FectoVIR®-AAV. Two independent experiments were conducted where each transfection was performed in duplicate. Three different plasmids ratio were evaluated (Transgene; Helper; RepCap). The % of full capsids showed is the results of qPCR:ELISA ratio (data provided by ABL).

