

# Quick and Easy Separation of Empty | Full AAV Particles Using Sartobind® Q Lab

Sartobind® Q Lab anion exchange membrane chromatography units can be used to separate empty from genome-filled adeno-associated virus (AAV) particles. Compared to classical resin columns, these membrane adsorber units allow for significantly higher flow rates to be used. In this case study, we summarize an empty | full separation process for AAV8 particles, although it should be noted that this can also be applied to other serotypes. Our step gradient elution method showed a 5.3-fold enrichment of full particles, making Sartobind® Q Lab an ideal choice for empty | full separation in AAV purification workflows.

## Materials

- Affinity-purified AAV sample
- Sartobind® Q Lab (0.41 mL membrane volume)
- Purification buffers (50 mM Tris-HCl pH 8.5 with 2 – 50 mM MgCl<sub>2</sub>)
- Liquid chromatography system or syringes
- AAV reference standards
- Octet® R8 BLI system with AAVX Biosensors
- dPCR system



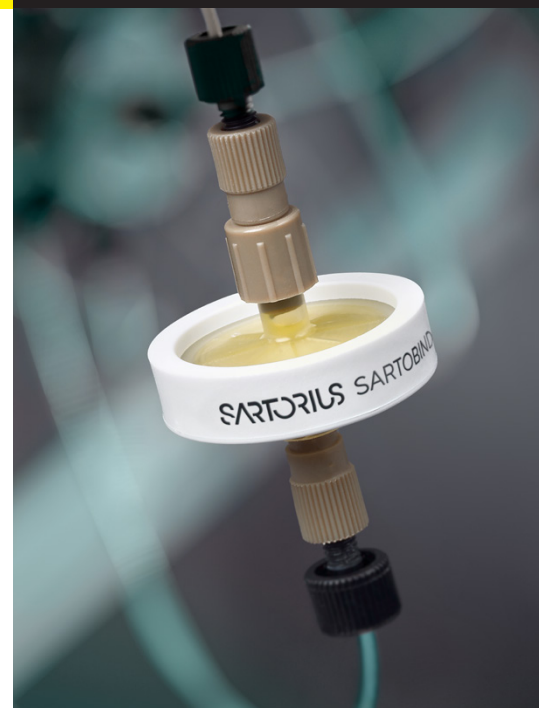
## Case Profile

### Objective:

This case study shows how Sartobind® Q Lab can be used to improve the efficiency of empty | full separation during the research and development of new AAV-based gene therapies. High flow rates and low shear forces speed up the polishing step used to enhance AAV purity.

### Keywords:

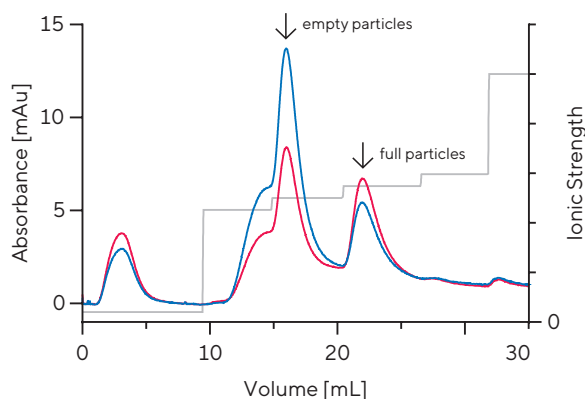
AAV, adeno associated virus, empty | full separation, full capsid enrichment, gene therapy, membrane chromatography, Octet®, Sartobind® Lab, anion exchange chromatography



# Results

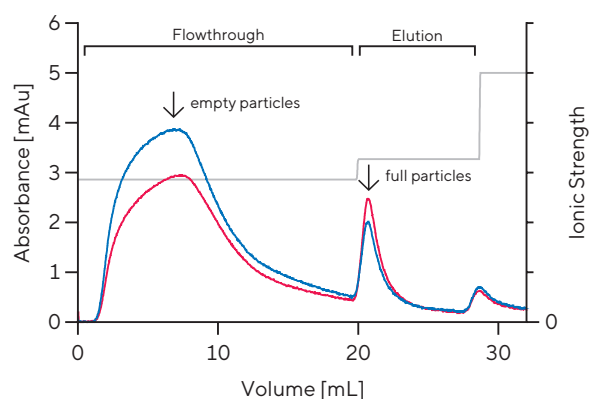
Ion exchange chromatography is commonly used to enrich genome-filled AAVs by exploiting the difference in isoelectric point between empty and full particles. **Figure 1** shows a 5.3-fold enrichment of full AAV8 particles when Sartobind® Q Lab was used in combination with a step elution gradient, where the loaded material and second elution peak contained 12.1% and 63.9% full particles, respectively.

## 1. Step Elution E | F Separation



To simplify the empty | full separation process, a flowthrough approach was developed. Sartobind® Q Lab was equilibrated at an ionic strength at which only full AAV8 particles would be captured during loading. **Figure 2** shows that loading and elution at 25 mM and 32.5 mM MgCl<sub>2</sub>, respectively, resulted in a single elution peak containing 44.9% full particles – a 3.7-fold enrichment factor.

## 2. Flowthrough E | F Separation



# Conclusion

Our results show that Sartobind® Q Lab can be effectively used for empty | full separation of AAV particles. Step elution and flowthrough approaches resulted in up to 5.3-fold enrichment of full AAV8 particles, relative to the affinity-purified starting material. Further data for AAV8 and other serotypes are available in our Application Note: [Establishing a Small Scale AAV Empty | Full Separation Process using Sartobind® Q Lab](#).

These methods can increase empty | full separation efficiency because the convective Sartobind® Q membrane supports faster flow rates than diffusive, resin-based columns. In addition, the flowthrough approach is easy to perform using syringes, eliminating the additional time and effort required to set up a liquid chromatography system. Therefore, Sartobind® Q Lab offers distinct advantages to accelerate and simplify empty | full separation of AAVs in research and development applications.

## Germany

Sartorius Lab Instruments GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Göttingen  
Phone +49 551 308 0

## USA

Sartorius Corporation  
3874 Research Park Drive  
Ann Arbor, MI 48108  
Phone +1 734 769 1600

 For further information, visit  
[sartorius.com](https://www.sartorius.com)