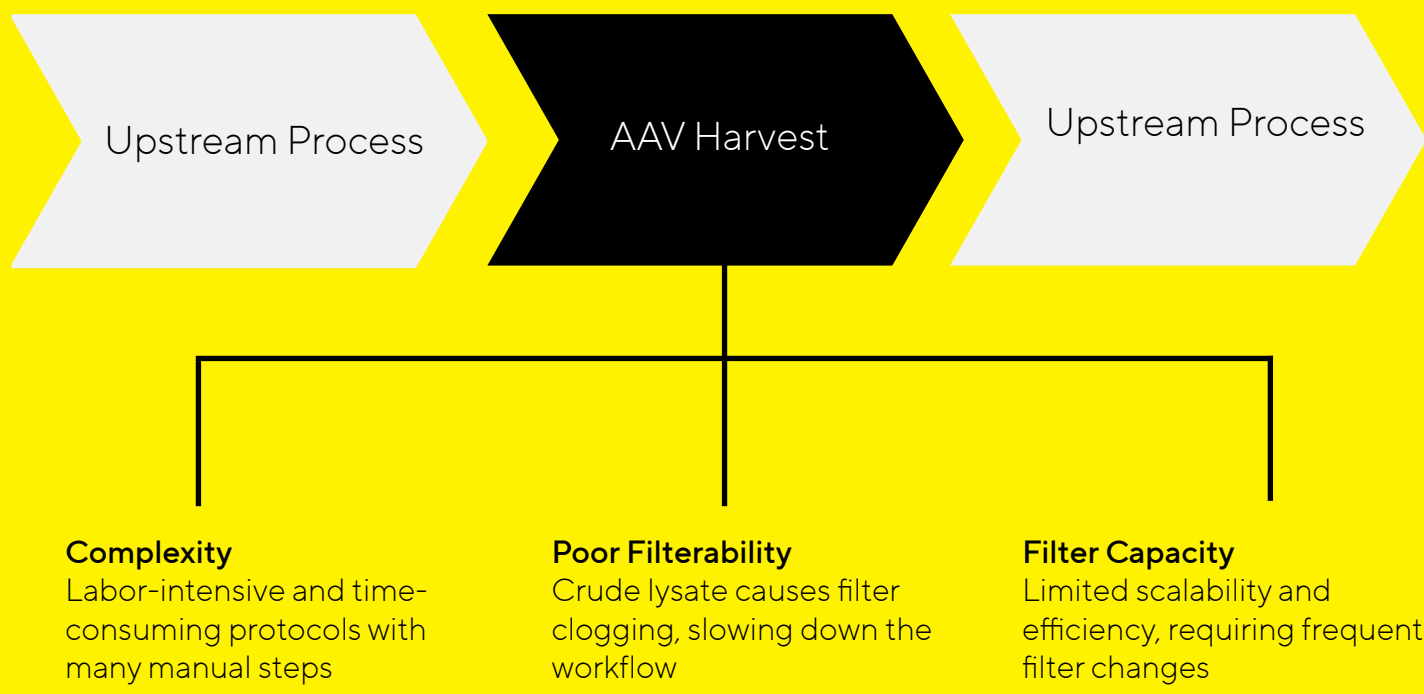


# Efficient Adeno-Associated Viral Vector (AAV) Clarification with Diatomaceous Earth

After production of AAV particles in host cells (e.g., HEK293 cells), you need to remove cell debris and other impurities from the crude lysate. This is necessary prior to purification and formulation of the AAV for therapeutic applications.

## AAV Harvest Pain Points



## Benefits of a Filter Aid

Filter aids such as diatomaceous earth (DE) improve filtration efficiency by preventing filter clogging. They are added to the filter medium (pre-coat) or to the fluid being filtered.



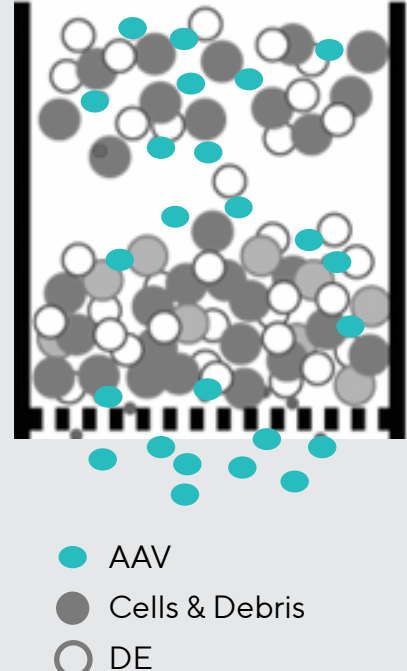
Diatomaceous Earth (DE)

### What is DE?

- Fossilized remains of diatoms
- Highly porous
- Insoluble & inert

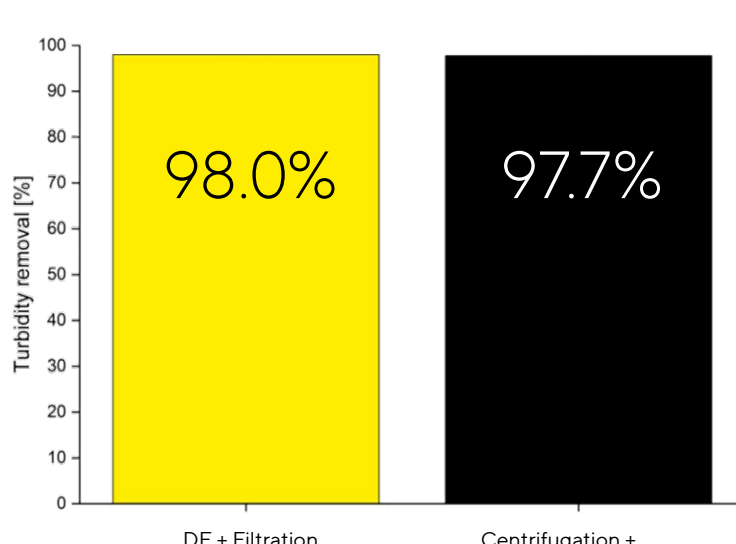
## How Does DE Filtration Work?

- DE grade C300 is designed for removal of mammalian cells and cell debris
- DE forms a porous layer that separates cells and debris
- AAV passes through easily
- One step clarification - no centrifugation

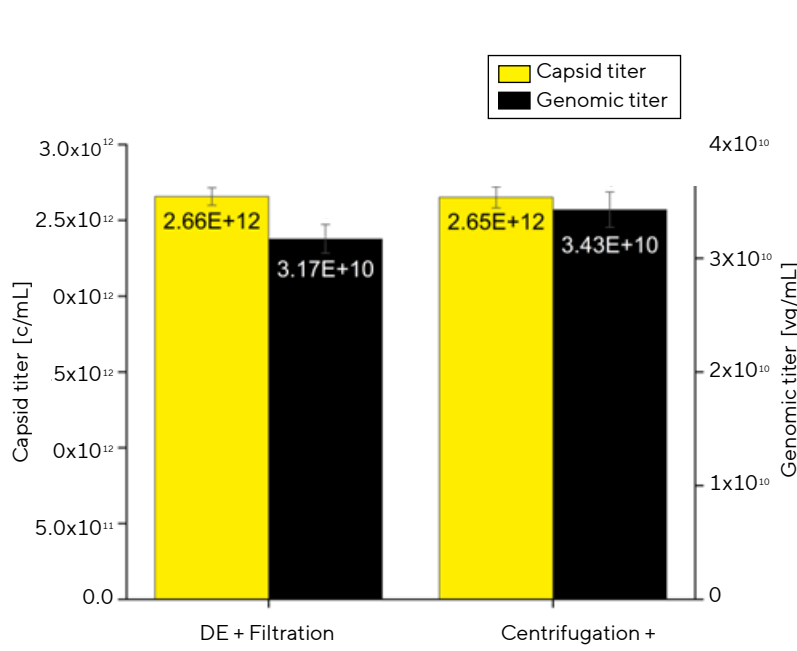


## One-Step AAV Harvest: Sartoclear Dynamics® Lab Filter Aid With DE

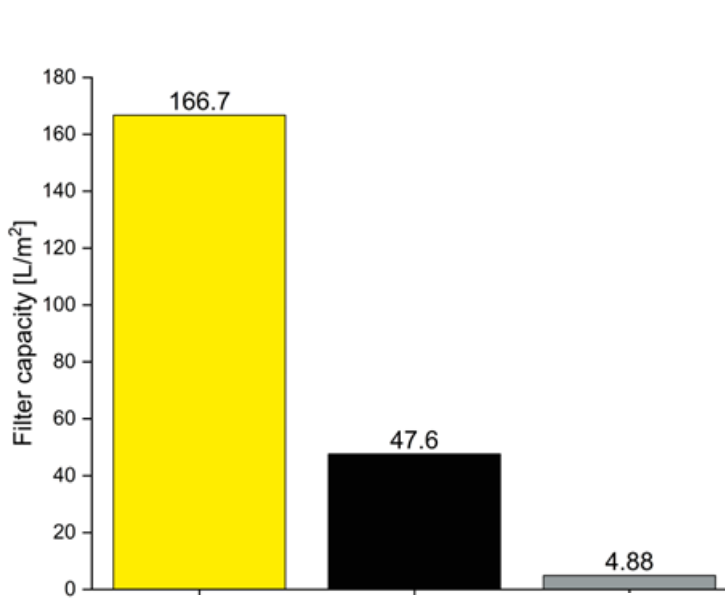
Same turbidity reduction as conventional methods



95% AAV5 (shown) & AAV8 recovery, (80% AAV2 recovery)



3.5-fold increase in filter capacity



Discover Sartoclear Dynamics® Lab Convenient Mammalian Cell Culture Harvesting