



Sartopore[®] platinum

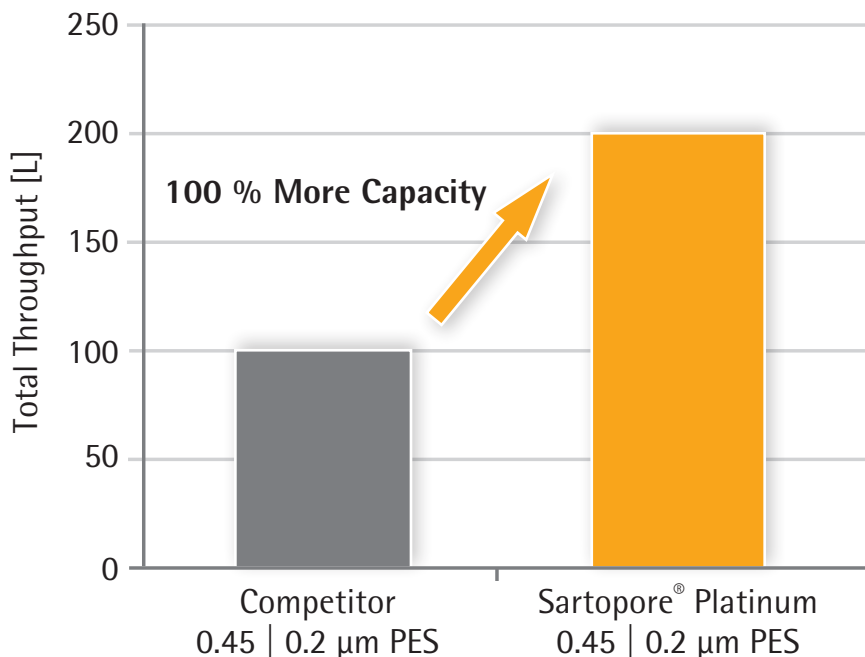
A New Vision of Perfection.
A New Class of Sterile Filtration.

Unique Performance in Filtration of Ophthalmics

Sartopore[®] Platinum is characterized by a unique and patented surface modification of the 0.45|0.2 µm rated Polyethersulfone (PES) membranes. In contrast to other PES filters in the market it is the highly hydrophilic low adsorptive membrane of Sartopore[®] Platinum which leads to an outstanding combination of high filtration capacity, high flow rate and low binding. This leads to lowest filter consumption and lowest adsorption of e.g. Benzalconiumchloride (BAC) and other preservatives. Sartopore[®] Platinum is the top-performer in the market for any Ophthalmic solutions.

Customer Example

- Application: ▶ Sterile filtration of a 0.5 % hypromellose solution used as an ophthalmic lubricant
- Challenge: ▶ Early blocking of existing filters and long filtration time due to high viscosity
- Customer's requirements: ▶ Reliable filtration despite high viscosity of hypromellose
▶ Reduce process times compared to current filter supplier
▶ Optimization of production costs
- Status with competition: ▶ 20" filter cartridge (PES, 0.45 | 0.2 µm) to filter 200 L in 2 hours
- Our solution: ▶ 10" Sartopore[®] Platinum to filter 200 L in less than 1 hour
- Customer benefits: ▶ >50 % reduction of process time
▶ 50 % reduction of filter consumption
▶ Costs savings of 150 US\$ per batch (75.000 US\$ per year)



Filtration Time

-50%

Filter Consumption

-50%