

The Proof

Powerful Filter Combination: Prefilter Sartoquard | Sterile Filter Sartopore® 2 XLM

Testfiltration at customer side lead to significant savings on pre and sterile filters.

Media Preparation

O.1 μm

Prefiltration

Final Filtration

Bioreactor

State with competition: 38×30 " prefilter with 0.1 μ m + 6×30 " final filter 0.1 μ m

Our solution: 12×30 " Sartoguard PES 0.1 μ m + 3×30 " Sartopore $^{\circ}$ 2 XLM (0.2 μ m + 0.1 μ m)

Customer benefit: Saving of **1.740** 30" filters per year (60 batches)

(together with other media steps even 2.580 30" filters per year)

