

## Adsorptive Pre-Filtration to Increase Virus Filter Performance and Overall Process Robustness in Blood Derived Processes

Volkmar Thom, Björn Hansmann, Jörg Hosch, Benjamin Schneider

Sartorius-Stedim Biotech GmbH, Göttingen, Germany, Membrane R&D

### Adsorptive pre-filters

The evaluation of virus filters is not confined only to their capacity to retain viruses. Indeed, selection of a virus filter is influenced by numerous factors. One factor gaining increase importance is process economics. Different adsorptive pre-filters have been introduced to the market for capacity increase of virus-retentive filters. Today's established adsorptive pre-filters are compared in the table below.

Depth Filter	CEX Membrane	Virosart® Max <sup>1</sup>
● Nearly independent of conductivity	● Affected by process conditions (pH, conductivity)	● Performance independent from process conditions (conductivity)
● High extractable   particle load	● Low extractable   particle load	● Low extractable   particle load
● Integrity test not available	● Integrity test not available	● Integrity test by air diffusion

<sup>1</sup> Sartorius patent DE102011105525-B4; US, EP and WO patents pending.  
<sup>2</sup> Method for removing biopolymer aggregates and viruses from a fluid<sup>2</sup>

### Characteristics of Virosart® Max

#### Working principle

- Combination of adsorptive capacity and size exclusion leads to removal of virus filter foulants
- Aggregates and/or small hydrophobic molecules are typical virus filter foulants

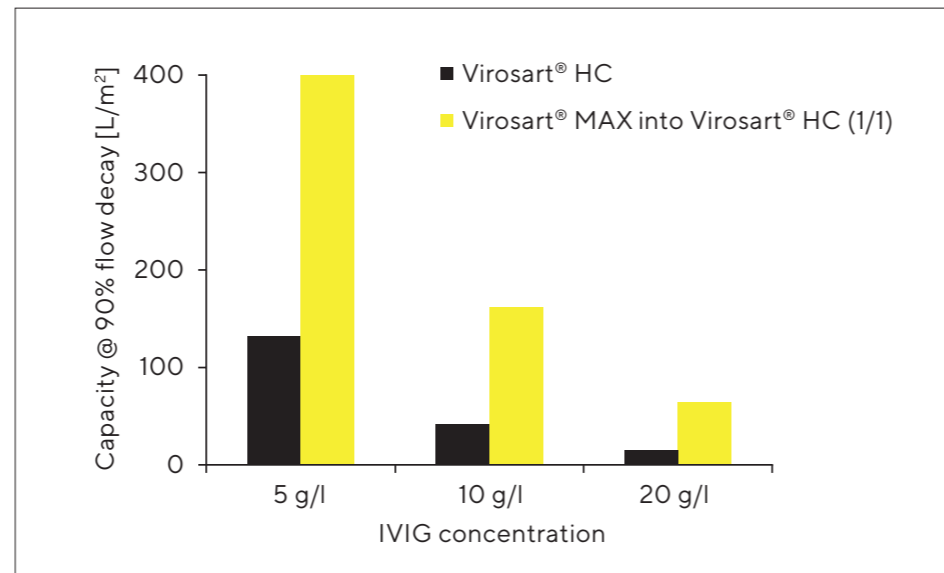
#### Filter Configuration

- Material: Optimized polyamide
- Pore size: 0.1 µm (nominal)
- Format: Triple-layer pleated elements
- Size: Available from 5 cm<sup>2</sup> to 30" elements



#### Higher capacity through aggregate reduction

The impact of Virosart® Max on the filtration of different IVIG concentrations (5, 10 and 20 g/L) through Virosart® HC 20 nm virus filter (5 cm<sup>2</sup> Minisart® devices) was analyzed. Filtrations have been performed with and without the use of Virosart® Max at 2.0 bar | 30 psi filtration pressure. Results were compared at 90% flow decay.

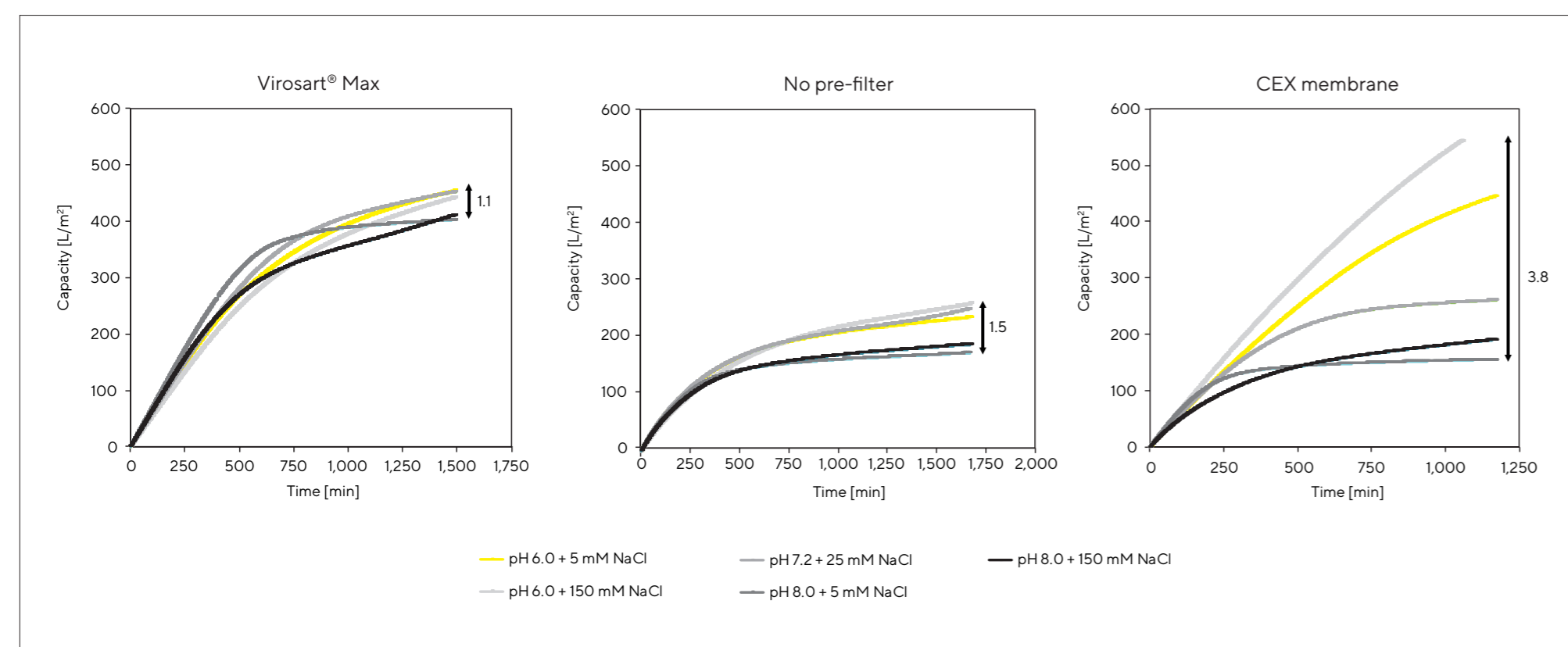


As a result, filtration capacity scales with solution concentration because the concentration of membrane fouling impurities scales accordingly.

#### Robust against process conditions

The effect of different pre-filtration strategies was evaluated for IVIG (5 g/L) in different buffer conditions at varying pH and ionic strength using Virosart® HC 20 nm virus filter (5 cm<sup>2</sup> Minisart® devices) at 2.0 bar | 30 psi.

As a result, the use of Virosart® Max results in lowest performance spread by varying process conditions.

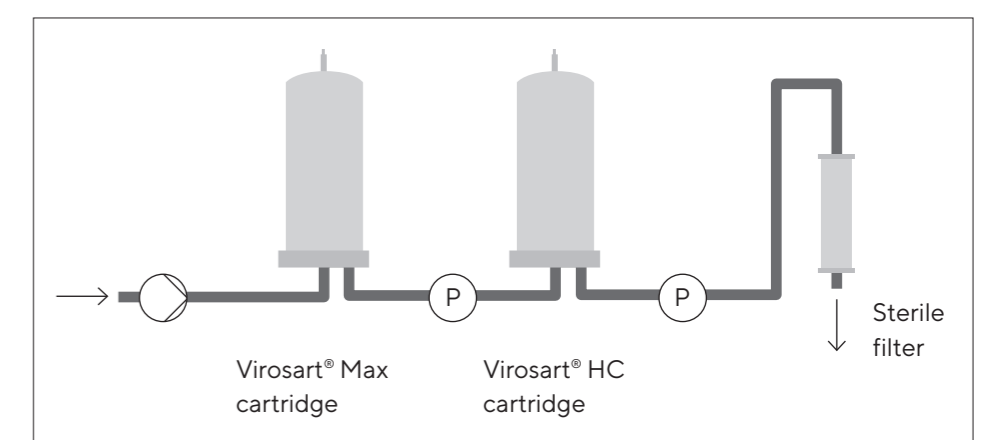


### Implementation

Cartridges and capule format of the filter allows flexible process implementation:

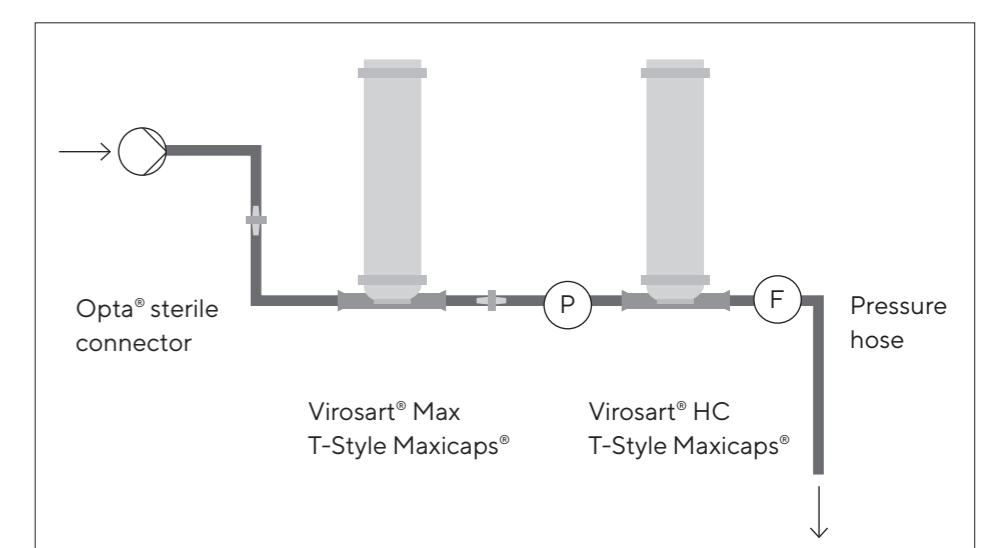
#### Stainless steel housing setup

- Robust setup
- Steam sterilization and pre-use integrity testing possible



#### Single-use setup

- Ease of use
- Flexible
- Pre-use integrity testing limited under fully-contained sterile conditions



#### Automated setup

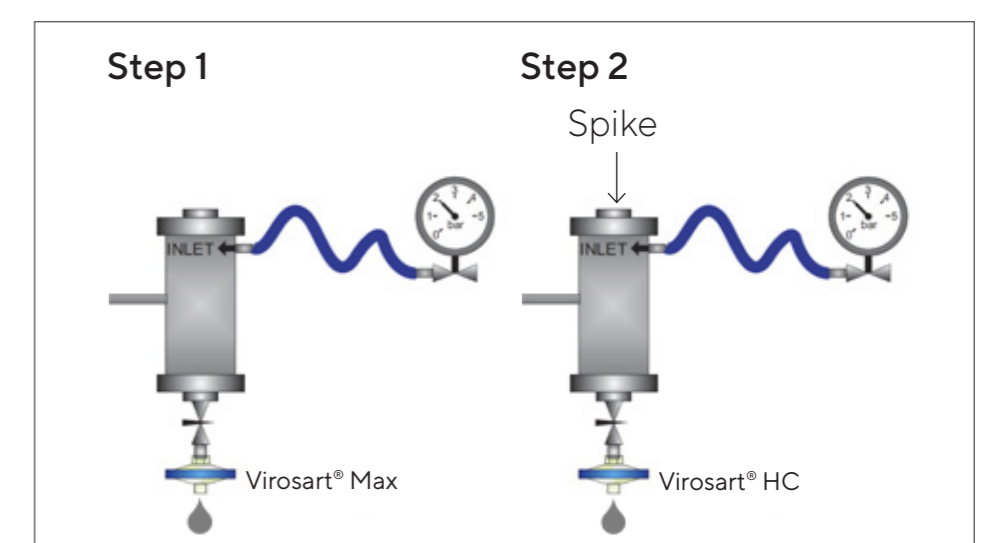
- Customized set-up
- High level of automation



### Spiking studies

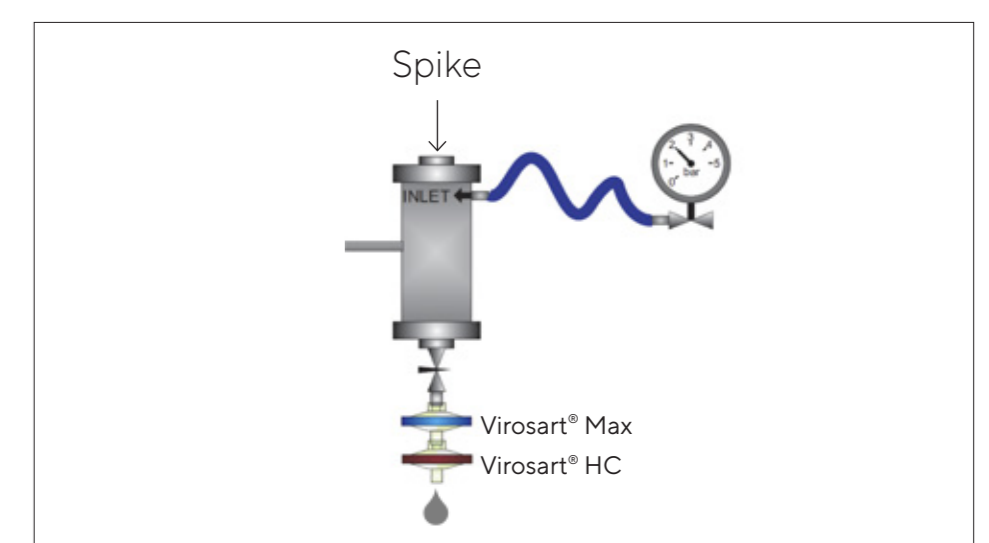
#### Preferred Option:

- Off-line pre-filtration (decoupled)
- Product is pre-filtered off-line and afterwards virus spike is added to the product feed
- Pressure | flow adaption over pre-filter
- Low capacity of virus filter by highly fouling feed streams
- Common approach in the industry
- Pre-filtration before validation to restore sample



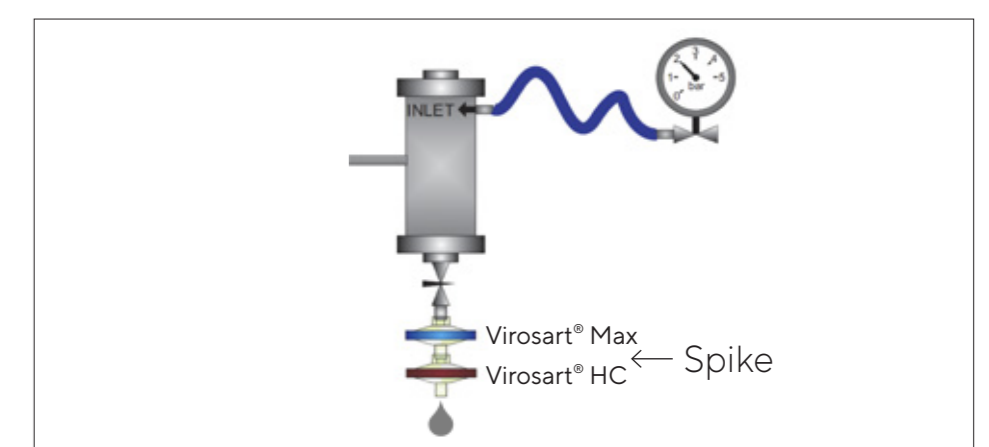
#### Alternative 1:

- In-line pre-filtration (coupled)
- Pre-filter and virus filter are run in-line and virus spike is added in-line.
- Virus retention by pre-filter not rated as robust
- Possible if pre-filter is tested independently for virus retention



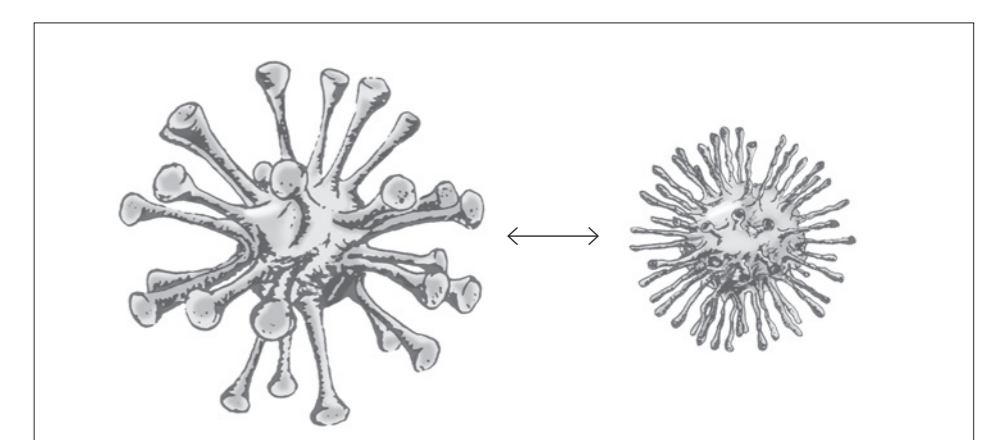
#### Alternative 2:

- In-line pre-filtration with in-line spiking
- Pre-filter and virus filter are run in-line, but the virus spike is added in-line after the pre-filter.
- Complex setup
- Difficult control of feed titer



#### Alternative 3:

- Spiking virus selection
- Validate virus-retentive filter for parvoviruses (PPV, MVM) and imply sufficient LRV for larger viruses (MuLV, PRV)
- Accepted by regulatory authorities?



### References

'Artifacts of Virus Filter Validation', P. Genest, H. Ruppach, C. Geyer, M. Asper, J. Parrella, B. Evans, A. Slocum, BioProcess International 2013.