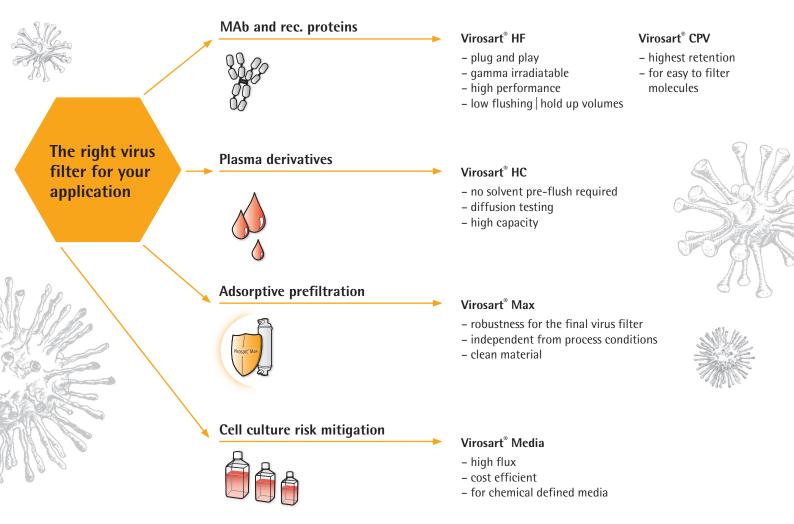


Virosart[®] virus filters Reliable retention under all circumstances



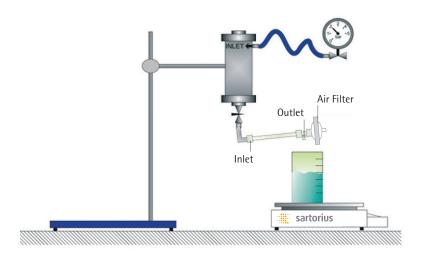
No impact on virus retention through pressure variations, high loads or process interruptions.

- ≥ 4 log for small non-enveloped viruses
- ≥ 6 log for large enveloped viruses



How to use Virosart® HF

Use Virosart® HF for virus filtration at the end of the purification process.





- install the device in the direction of filtration
- connect vent filter in a slightly upward position (see picture)

Wetting

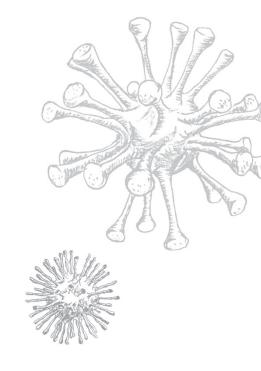
- rinse filter with buffer | WFI in the direction of filtration at an inlet pressure of 2 bar for 15 min
- typical flow rates with WFI:
 200 LMH ± 25% @ 1 bar, 25°C

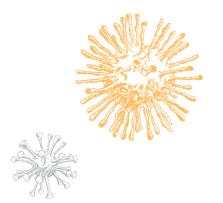
3 Filtration

- replace buffer | WFI by the solution to be filtered
- keep vent filter connected during the filtration
- start filtration at the desired filtration pressure

Max. allowable differential pressure

- direction of filtration: at 20°C max. 3 bar | 43.5 psi (up to 5 bar with plug instead of air filter)
- reversed direction of filtration: at 20°C max. 2.5 bar | 36.3 psi







For detailed description please visit: www.sartorius-stedim.com/virus-filtration or scan QR code.