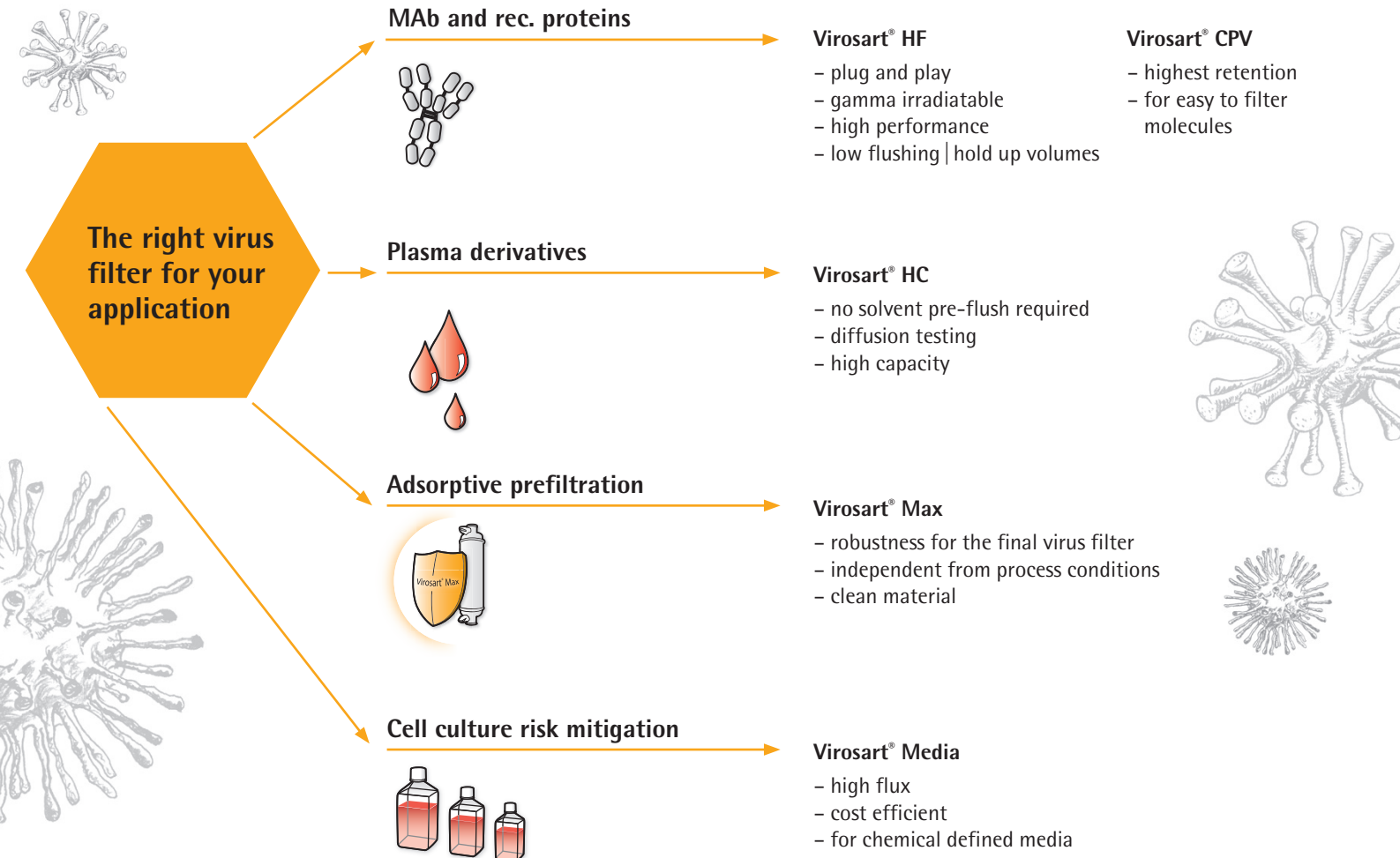


# Virosart® virus filters

## Reliable retention under all circumstances

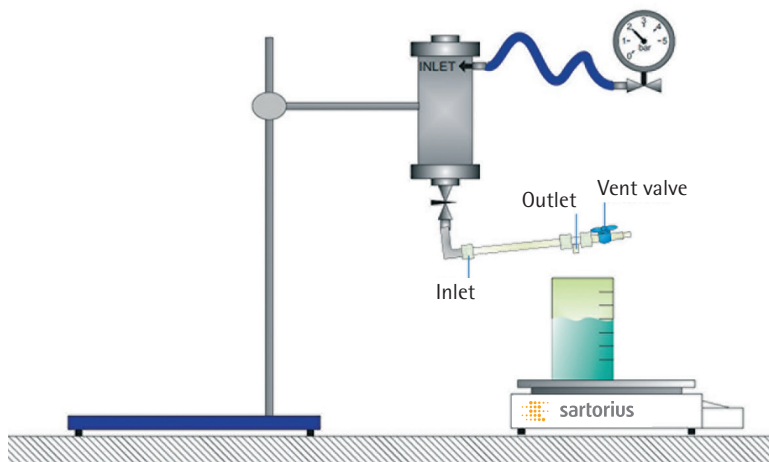


**No impact on virus retention through pressure variations, high loads or process interruptions.**  
 $\geq 4$  log for small non-enveloped viruses  
 $\geq 6$  log for large enveloped viruses



# How to use Virosart® Media

Use Virosart® Media for virus filtration upstream of the bioreactor for chemically defined cell culture media.



## 1 Installation

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

## 2 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:  
450 LMH  $\pm$  25% @ 1 bar, 25°C

## 3 Filtration

- replace buffer | WFI by the solution to be filtered
- open the vent valve after every change within the reservoir
- start filtration at the desired filtration pressure

### Max. allowable differential pressure

- direction of filtration: at 20°C max. 5 bar | 73 psi
- reversed direction of filtration:  
at 20°C max. 1 bar | 14.6 psi



For detailed description please visit:  
[www.sartorius-stedim.com/virus-filtration](http://www.sartorius-stedim.com/virus-filtration)  
or scan QR code.

