

## Virosart® Virus Filters

Robust, Efficient and Reliable

State-of-the-art virus retention, ensuring performance even under challenging process conditions such as pressure fluctuations, high protein loads, and process interruptions.

$\geq 4 \log_{10}$  for small non-enveloped viruses

$\geq 6 \log_{10}$  for large enveloped viruses



# The Right Virus Filter for Your Application

| Application                   | Downstream                                                                                                                                                                                                                                                                                                                                                               |
|-------------------------------|--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| mAbs and recombinant proteins | <div><p><b>Virosart® HF</b></p><ul style="list-style-type: none"><li>▪ High flux</li><li>▪ Sterile delivery</li><li>▪ Plug and play</li><li>▪ Low flushing   hold up volumes</li></ul><p><b>Virosart® CPV</b></p><ul style="list-style-type: none"><li>▪ Caustic stable</li></ul></div> |
| Plasma derivatives            | <div><p><b>Virosart® HC</b></p><ul style="list-style-type: none"><li>▪ High capacity</li><li>▪ Robustness   low blocking</li></ul></div>                                                                                                                                                |
| Prefiltration                 | <div><p><b>Virosart® Max</b></p><ul style="list-style-type: none"><li>▪ Adsorptive   size exclusion</li><li>▪ Aggregate removal</li><li>▪ Protect final virus filter</li></ul></div>                                                                                                    |
| Application                   | Upstream                                                                                                                                                                                                                                                                                                                                                                 |
| Cell culture media            | <div><p><b>Virosart® Media</b></p><ul style="list-style-type: none"><li>▪ Risk Mitigation</li><li>▪ High flux</li><li>▪ For chemically defined media</li></ul></div>                                                                                                                  |



## Find out more

Discover how Virosart® can enhance  
your filtration process:  
[sartorius.com/virus-filtration](https://sartorius.com/virus-filtration)

## Germany

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen  
Phone +49 551 308 0

