

# SARTORIUS

## Virosart® Virus Filters Reliable Retention Under All Circumstances

**No impact of pressure variations,  
high loads and process interrup-  
tions on virus retention.**

≥ 4 log for small non-enveloped  
viruses

≥ 6 log for large enveloped viruses



# The Right Virus Filter for Your Application

---

## Application

## Virosart® Virus Filter

MAb and rec. proteins



Virosart® HF

- plug and play
- gamma irradiatable
- high performance
- low flushing | hold up volumes



Virosart® CPV

- highest retention
- for easy to filter molecules

---

Plasma derivatives



Virosart® HC

- no solvent pre-flush required
- diffusion testing
- high capacity

---

Adsorptive prefiltration



Virosart® Max

- robustness for the final virus filter
- independent from process conditions
- clean material

---

Cell culture risk mitigation



Virosart® Media

- high flux
- cost efficient
- for chemically defined media

---

## Find out more

For detailed description please visit:  
[www.sartorius.com/virus-filtration](http://www.sartorius.com/virus-filtration)

