

Virosart[®] Virus Filters Reliable Retention Under Challenging Conditions

**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

- High 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 information please visit:
www.sartorius.com/virus-filtration

Germany

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

