



**sartorius stedim**  
biotech

# Highest virus safety for your patient

## Next generation virus clearance concepts



**The right solution for every application**

**Virosart®**

The 20 nm PES filters for robust and efficient virus removal

**Sartobind®**

Membrane chromatography for virus & contaminant clearance



[www.sartorius-stedim.com/virus-clearance](http://www.sartorius-stedim.com/virus-clearance)



## 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



**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**

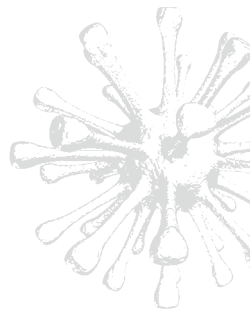
- robustness for the final virus filter
- independent from process conditions
- clean material

**Cell culture risk mitigation**



**Virosart® Media**

- high flux
- cost efficient
- for chemical defined media



Learn more about our Virosart® virus filters at [www.sartorius-stedim.com/virus-filtration](http://www.sartorius-stedim.com/virus-filtration) or scan QR code.