

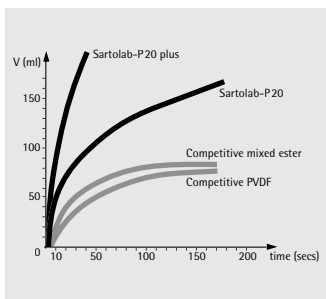
Sartolab® P20 and Sartolab® P20 plus for Reliable Sterile Filtration of Tissue Culture Solutions



Ready-to-use units which, attached to a membrane pump or tube pump, quickly and reliably sterilize 100 ml to 5 liter of media and aqueous solutions.

The combination of a large filtration area (20 cm²) and an automatic de-aeration ensures high flow rates and optimal total throughputs. Automatic venting of any trapped air through the PTFE membrane-protected vent ports ensures that the entire filter surface is used for effective filtration.

Sartolab P20 units have an extraordinarily favourable price-performance ratio. Often, the total filterable volume can even be doubled due to an integrated binder-free glass fibre prefilter.



Top part: results with contaminated medium (DMEM + 10% FCS), at 1 bar differential pressure.

Specifications for Sartolab P20 units

Connectors	Inlet, luer lock inner cone or 6–12 mm stepped hose nipple. Outlet, 6–12 mm stepped hose nipple
Biosafety	Pass the USP Plastics-Class VI-Test
Bubble point	Min. value with water, 3.2 bar (320 kPa, 46 psi)
Flow rate	For water 250 ml/min at $\Delta p = 1$ bar (100 kPa, 14.5 psi)
Filtration area	20 cm ²
Filling volume	6 ml
Housing burst pressure	> 5 bar (500 kPa, 72.5 psi)
Materials	Cellulose acetate membrane filter (0.2 μ m). PTFE airfilter. Polycarbonate housing
Max. recommended inlet pressure	3 bar (300 kPa, 43.5 psi)
Protein adsorption	Less than 10 μ g γ -Globulin/cm ²
Hold-up volume	0.3 ml after (1.3 ml before) bubble point
Toxicity	Non-toxic as confirmed with L929 fibroblast cells of mice and with MRC-5 lung cells of human embryonic origin
Accessories	Integrity holder 18099

Specifications for Sartolab P20 plus units

As for P20, except	
Filling volume	5.5 ml
Materials	Supplemented with a binder-free glass fibre prefilter
Protein adsorption	Varies due to the prefilter
Hold-up volume	0.9 ml after (1.8 ml before) bubble point

Order numbers see next page.

**Order numbers for Sartolab P20 units**

18052 D	With hose nipple inlet connection, pack of 10
18053 D	With luer lock inlet connection, pack of 10

**Order numbers for Sartolab P20 plus units**

18056 D	With hose nipple inlet connection, pack of 10
18058 D	With luer lock inlet connection, pack of 10

Recommended accessories are described on page 72.
Special brochure available on request. Order no. SL-3009-e