

Sartolab® P20 Portfolio Expansion

All new part numbers are highlighted in yellow.

Sartolab® P20

Order Number	Filter Material	Inlet	Outlet	Sterilization (EO)	Qty./Pack
18064-----D	0.2 µm PES	Female Luer-Lock	Male Luer-Lock	yes	10
18075-----D	0.2 µm PES	Female Luer-Lock	Hose barb	yes	10
18075-----UPN	0.2 µm PES	Female Luer-Lock	Hose barb	no	100
18089-----D	0.2 µm PES	Hose barb	Hose barb	yes	10
18090-----D	0.2 µm PES	Female Luer-Lock	Male Luer-Lock	yes	10

Sartolab® P20 Plus

Order Number	Filter Material	Inlet	Outlet	Sterilization (EO)	Qty./Pack
18068-----D	Quartz microfibers & 0.2 µm PES	Female Luer-Lock	Hose barb	yes	10
18076-----N	Quartz microfibers & 0.45 µm PES	Hose barb	Hose barb	no	100
18091-----D	Quartz microfibers & 0.2 µm PES	Hose barb	Hose barb	yes	10
18092-----D	Quartz microfibers & 0.2 µm PES	Female Luer-Lock	Male Luer-Lock	yes	10

Sartolab® P20 Prefilter

Order Number	Filter Material	Inlet	Outlet	Sterilization (EO)	Qty./Pack
18072-----D	Quartz microfibers	Female Luer-Lock	Hose barb	no	10

 New part numbers