



미생물 품질시험의 기준

'Sartorius Vacuum Filtration System'

소모품 구매 시

300만원 상당의 3 Way Microsart® Manifold

무상 제공

*소모품 : Microsart ADD filter (20ea)

*기간 : 4월 22일 - 9월 30일

신청하기 ▶



Microsart® Manifold

FREE!



- 미생물 분석, 무균시험, 및 미세 잔류물 분석 등 다양한 Application 활용 가능
- 다양한 Filter Holder 장착 가능
- 교차 오염을 최소화 시키는 Venting Parts와 Autoclave 가능한 Full SS 316L 재질
- 낮은 높이(9.8cm)로 실험 편의성 증가



Ordering Information

Microsart® @Filter With CN Filter

Microsart® @Filter 100, with lid, packaged on trays, ideal for use in clean benches

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.2 µm	CN white black	100 mL	24	16D01-10-07-TG
0.45 µm, High Flow	CN white black	100 mL	24	16D01-10-H6-TG
0.45 µm, High Flow	CN gray white**	100 mL	24	16D03-10-H6-TG
0.45 µm	CN green dark green	100 mL	24	16D02-10-06-TG

Microsart® @Filter 250, with lid, packaged on trays, ideal for use in clean benches

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.2 µm	CN white black	250 mL	16	16D01-25-07-TF
0.45 µm, High Flow	CN white black	250 mL	16	16D01-25-H6-TF
0.45 µm, High Flow	CN gray white**	250 mL	16	16D03-25-H6-TF
0.45 µm	CN green dark green	250 mL	16	16D02-25-06-TF
0.65 µm	CN gray white**	250 mL	16	16D03-25-05-TF

Microsart® @Filter 100, stacked and packaged in bags, ideal for use with Microsart® Funnel Dispenser

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.2 µm	CN white black	100 mL	60	16D01-10-07-BL
0.45 µm, High Flow	CN white black	100 mL	60	16D01-10-H6-BL
0.45 µm, High Flow	CN gray white**	100 mL	60	16D03-10-H6-BL
0.45 µm	CN green dark green	100 mL	60	16D02-10-06-BL

Microsart® @Filter 250, stacked and packaged in bags, ideal for use with Microsart® Funnel Dispenser

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.2 µm	CN white black	250 mL	48	16D01-25-07-BK
0.45 µm, High Flow	CN white black	250 mL	48	16D01-25-H6-BK
0.45 µm, High Flow	CN gray white**	250 mL	48	16D03-25-H6-BK
0.45 µm	CN green dark green	250 mL	48	16D02-25-06-BK
0.65 µm	CN gray white**	250 mL	48	16D03-25-05-BK

Microsart® @Filter 100, individually sterile packaged, with lid

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.45 µm, High Flow	CN white black	100 mL	27	16D01-10-H6-ACG
0.45 µm, High Flow	CN black white	100 mL	27	16D03-10-H6-ACG
0.2 µm	CN white black	100 mL	27	16D01-10-07-ACG

Microsart® @Filter 250, individually sterile packaged, with lid

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.45 µm, High Flow	CN white black	250 mL	18	16D01-25-H6-ACF
0.45 µm, High Flow	CN black white	250 mL	18	16D03-25-H6-ACF
0.2 µm	CN white black	250 mL	18	16D01-25-07-ACF

Microsart® @Filter With PVDF Filter

Microsart® @Filter 100, stacked and packaged in bags, ideal for use with Microsart® Funnel Dispenser

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.45 µm	PVDF white no grid	100 mL	60	16D04-10-06-BL***

Microsart® @Filter 100, individually sterile packaged, with lid

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.45 µm	PVDF white no grid	100 mL	27	16D04-10-06-ACG***

Microsart® @Filter 100, with lid, packaged on trays, ideal for use in clean benches

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.45 µm	PVDF white no grid	100 mL	24	16D04-10-06-TG***

Microsart® @Filter With Polycarbonate Filter

Microsart® @Filter 100, stacked and packaged in bags, ideal for use with Microsart® Funnel Dispenser

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.45 µm	PC white no grid	100 mL	60	16D06-10-06-BL***

Microsart® @Filter With Regenerated Cellulose Filter

Microsart® @Filter 100, with lid, packaged on trays, ideal for use in clean benches

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.45 µm	RC white (w/o grid)	100 mL	24	16D05-10-06-TG***

Microsart® @Filter 100, stacked and packaged in bags, ideal for use with Microsart® Funnel Dispenser

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.45 µm	RC white (w/o grid)	100 mL	60	16D05-10-06-BL***



Find out more

사토리우스코리아바이오텍(유)
경기도 성남시 분당구 판교역로 220 플리드스페이스 8층
E. marketing@sartorius.co.kr
www.sartorius.co.kr