

# SARTORIUS

많은 관심과 성원에 힘입어 진행하는

## Bottle Top Filter 앙코르프로모션!

Multistation 무상 증정!



바틀탑필터 12박스 + Storage bottle 12박스 구매 시,  
Hands Free 필터링 시스템 무상 증정

### Your Key Benefits

- 무상 증정 (약 270만원 상당)  
50, 150, 250, 500, 1000ml 필터 호환  
최대 6개까지 동시에 여과할 수 있는  
Hands Free 바틀탑필터 거치대
- 프로모션 대상 제품

품명	품번
180E04-E	Sartolab RF 500, 0.22um
180E05-E	Sartolab RF 1000, 0.22um
180F04-E	Sartolab RF 500, 0.45um
180F05-E	Sartolab RF 1000, 0.45um
180-22-E	Sartolab Bottle 150ml
180-23-E	Sartolab Bottle 250ml
180-24-E	Sartolab Bottle 500ml
180-25-E	Sartolab Bottle 1000ml

기간 : 2023.12.31 까지  
(제품 소진 시 조기 마감)

※ 10대 한정수량

QR  
CODE

Find out more



Bottle top filter 500ml & 1L



Storage bottle  
150, 250, 500, 1000ml

## SARTORIUS의 Bottle top filter 제품

제품 종류	품명	품번
	180E01-2	Sartolab RF 50; PES; 0.22 µm; 24 pcs
	180E01-E8	Sartolab RF 50; PES; 0.22 µm; 96 pcs
	180E02-E	Sartolab RF 150; PES; 0.22 µm; 12 pcs
	180E03-E	Sartolab RF 250; PES; 0.22 µm; 12 pcs
	180E04-E	Sartolab RF 500; PES; 0.22 µm; 12 pcs
	180E05-E	Sartolab RF 1000; PES; 0.22 µm; 12 pcs
	180D03-E	Sartolab RF 250; PES; 0.1 µm; 12 pcs
	180D05-E	Sartolab RF 1000; PES; 0.1 µm; 12 pcs
	180F01-2	Sartolab RF 50; PES; 0.45 µm; 24 pcs
	180F01-E8	Sartolab RF 50; PES; 0.45 µm; 96 pcs
	180F02-E	Sartolab RF 150; PES; 0.45 µm; 12 pcs
	180F03-E	Sartolab RF 250; PES; 0.45 µm; 12 pcs
	180F04-E	Sartolab RF 500; PES; 0.45 µm; 12 pcs
	180F05-E	Sartolab RF 1000; PES; 0.45 µm; 12 pcs
		180E12-E
180E13-E		Sartolab BT 250; PES; 0.22 µm; 12 pcs
180E14-E		Sartolab BT 500; PES; 0.22 µm; 12 pcs
180E15-E		Sartolab BT 1000; PES; 0.22 µm; 12 pcs
180F15-E		Sartolab BT 1000; PES; 0.45 µm; 12 pcs
		180-22-E
	180-23-E	Sartolab bottle, 250 mL
	180-24-E	Sartolab bottle, 500 mL
	180-25-E	Sartolab bottle, 1000 mL

명함