

SARTORIUS

Bottle Top Filter 구매하고 추가혜택 받아보세요!



기간 : 2025. 09. 01 ~ 2025. 10. 31

*EVENT 1, 2 중 택 1



▲ 프로모션 신청하기

EVENT 1

Sartolab RF 5 박스 구매 시,
Multistation 30% 할인

(Sartolab RF : 150ml, 250ml, 500ml, 1000ml)



EVENT 2

Sartolab RF 1박스 구매 시,
Bottle 1박스 무상 제공

(동일 볼륨에 한함, 재고 소진 시 까지)



X 12
FREE