

Sartorius 수동피펫 무상증정 프로모션



싸토리우스 Tacta 프로모션 이벤트!

기간 : 2023.04.10 ~ 2023.09.30 까지



+



+



Your Key Benefits

- 수동피펫 싱글채널 3자루 구매시,
+ 싱글채널 1자루 무상증정
+ 스탠드 또는 Tip 1 box 선택 증정
(Tip 선택할 경우 5ml : 5 tray, 10ml : 2 tray)
- 수동피펫 멀티채널
15% 특별할인!!!



가까운 싸토리우스 대리점 또는 아래 연락처로 문의 주세요!

T. 031.622.5785

E. marketing@sartorius.co.kr

명함



Find out more

Tacta Mechanical Pipette



Perfectly Balanced



Channels	Volume Range (µl)	Order no.
1	0.1-3	LH-729010
	0.5-10	LH-729020
	2-20	LH-729030
	10-100	LH-729050
	20-200	LH-729060
	100-1,000	LH-729070
	500-5,000	LH-729080
8	1,000-10,000	LH-729090
	0.5-10	LH-729120
	5-100	LH-729130
12	30-300	LH-729140
	0.5-10	LH-729220
	5-100	LH-729230
	30-300	LH-729240

Tacta가 선사하는 4가지 Key points!



① 인체공학적 디자인과 가벼운 Pipetting Force



② 전체 고압증기멸균 가능 내구성 높은 재질



③ 정확한 볼륨 조절과 Locking 기능



④ Warranty 2년 제공



Find out more

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