

## Picus<sup>®</sup>2 사전예약 이벤트



신제품 런칭 기념 충전 스탠드 무상 증정!

기간: 2023.12.31 까지



1자루 구매

FREE!



3자루 구매

FREE!

\*30 set 한정

### Your Key Benefits

- 1자루 구매 시  
+ 1구 충전 스탠드 증정 (30만원 상당)
- 3자루 이상 구매 시  
+ 4구 충전 스탠드 증정 (80만원 상당)

### Picus 2 4 Key Point !!



Mobile app을 통한  
편리한 설정 변경



Application에 따른  
8가지 모드



100g의 가벼운 무게



1 시간 충전 시  
1,000 cycle 가능



Find out more

## Picus®2 Electronic Pipette

The Next Generation of Pipetting

Channels	Volume Range ( $\mu$ l)	Order no.
1	0.2-10	LH-747021
	5-120	LH-747041
	10-300	LH-747061
	50-1,000	LH-747081
	100-5,000	LH-747101
	500-10,000	LH-747111
8	0.2-10	LH-747321
	5-120	LH-747341
	10-300	LH-747361
	50-1,200	LH-747391
12	0.2-10	LH-747421
	5-120	LH-747441
	10-300	LH-747461
	50-1,200	LH-747491



가까운 사토리우스 대리점 또는 아래 연락처로 문의 주세요!

T. 031.622.5785

E. [marketing@sartorius.co.kr](mailto:marketing@sartorius.co.kr)

명함



Find out more