

SARTORIUS



새로운 도전과 시작 
싸토리우스가 응원 합니다!

Pipette
프로모션

~ 2024년 4월 30일 까지, 선착순 100 set 소진 시 종료

1 Basic Package

Tacta 3(자루) : (10,100,1000ul 또는 20,200,1000ul)

~~1,300,000 원~~ →

750,000 원

2 Advanced Package

Basic Package + Picus2 1(자루)

~~2,500,000 원~~ →

1,200,000 원

3 All-in-One Package

Advanced Package + Midi Plus

~~3,200,000 원~~ →

1,500,000 원

All-in-One Package

Advanced Package

Basic Package



Tacta 3



Picus2



Midi Plus(Pipette aid)



Free Gifts!!!

Tip 1 Box (10 tray) 무상증정!



10ul / 200ul / 1000ul 중 택 1

Picus 2 multi channel 변경 안내

- 25만원 추가 시 8-ch Multi Pipette 변경 가능
- 30만원 추가 시 12-ch Multi Pipette 변경 가능

Tacta[®] Mechanical Pipette

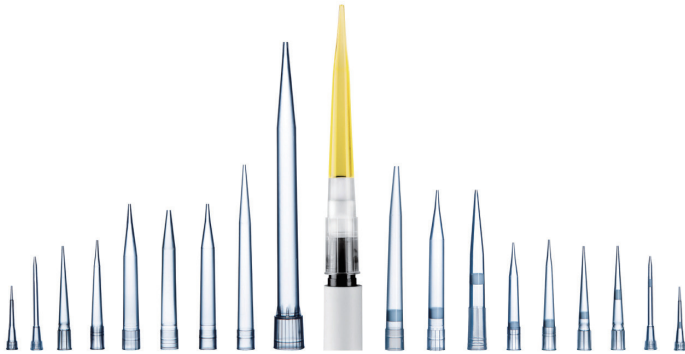


- 가볍고 부드러운 pipetting
- Volume lock 가능
- Autoclave 가능
- 다양한 Brand Tip 호환 가능

Picus2 Electronic Pipette



- 인체공학적인 디자인
- 다양한 피펫팅 모드
- 블루투스를 통한 앱 연동
- 샘플 손실 최소화



이번 프로모션을 통해 사토리우스의 고품질 팁을 무료로 경험해보세요!

Optifit Tips

- 샘플 손실을 최소화하는 고품질 팁
- 다양한 volume 팁 보유
- 다양한 Brand 피펫 호환 가능
- DNase, RNase, Endotoxin Free

궁금하신 사항은 아래 연락처로 문의 주세요!



Find out more

사토리우스코리아바이오텍(유)
 경기도 성남시 분당구 판교역로 220 쉐리드스페이스 8층
 E . marketing@sartorius.co.kr T.031.622.5700
 www.sartorius.co.kr

명함

선택 가능 모델

Order no.	Tacta (수동 Pipette) Pack 중 택 1
LH-729670	Tacta Pipette 3-pack 10(1ch-10/100/1000)
LH-729671	Tacta Pipette 3-pack 20(1ch-20/200/1000)

Order no.	Pipette Tip 무료 제공 제품 중 택 1
790011	Tip 10, Single Tray (10x96), Prs
790201	Tip 200, Single Tray (10x96), Prs,
791001	Tip 1000, Single Tray (10x96), Prs,



▪ Tacta 3 pack 구성품 : 피펫,스탠드,볼륨별 팁 1tray

Order no.	Picus2(전동 Pipette) 모델 중 택 1
LH-747021	Picus 2 Electronic Pipette, 1-ch, 0.2 - 10 ul
LH-747041	Picus 2 Electronic Pipette, 1-ch, 5-120 ul
LH-747061	Picus 2 Electronic Pipette, 1-ch, 10-300 ul
LH-747081	Picus 2 Electronic Pipette, 1-ch, 50-1000 ul
LH-747101	Picus 2 Electronic Pipette, 1-ch, 100-5000 ul
LH-747111	Picus 2 Electronic Pipette, 1-ch, 500-10000 ul

Order no.	Picus2(전동 Pipette) Multi Channel 변경 가능 품목
LH-747321	Picus 2 Electronic Pipette, 8-ch, 0.2-10 ul
LH-747341	Picus 2 Electronic Pipette, 8-ch, 5-120 ul
LH-747361	Picus 2 Electronic Pipette, 8-ch, 10-300 ul
LH-747391	Picus 2 Electronic Pipette, 8-ch, 50-1200 ul
LH-747421	Picus 2 Electronic Pipette, 12-ch, 0.2-10 ul
LH-747441	Picus 2 Electronic Pipette, 12-ch, 5-120 ul
LH-747461	Picus 2 Electronic Pipette, 12-ch, 10-300 ul
LH-747491	Picus 2 Electronic Pipette, 12-ch, 50-1200 ul

Order no.	Midi Plus (Pipette aid)
710931	Midi Plus, 1-100 ml with AC-adaptor



궁금하신 사항은 아래 연락처로 문의 주세요!

T. 031.622.5700

E. marketing@sartorius.co.kr

명함



Find out more