

## Picus®2 런칭 기념 피펫 프로모션



수동, 전동 상관 없이 무한 혜택 제공!

기간: 2023.12.31 까지

### EVENT 1 피펫 Trade in 프로모션

- 헌 피펫 반납하고 새 피펫 받아주세요!  
가지고 있는 피펫 반납 시(브랜드 무관)  
신규 피펫 30% 할인



Tacta



Picus 2



### EVENT 2 Tacta multi pack 할인

- 피펫 세트에 Stand와 Tip까지 함께  
알뜰하게 가져가세요!  
Tacta multi pack 제품 20% 할인  
(구성품은 하기 이미지 참조)



QR  
CODE

Find out more

# Sartorius Pipette

Perfectly Balanced

Event 1.  피펫 Trade in 프로모션	Channels	Volume Range ( $\mu\ell$ )	Tacta	Picus 2	
	1	3		LH-729010	-
		10		LH-729020	LH-747021
		20		LH-729030	-
		100		LH-729050	-
		120		-	LH-747041
		200		LH-729060	-
		300		-	LH-747061
		1,000		LH-729070	LH-747081
		5,000		LH-729080	LH-747101
10,000			LH-729090	LH-747111	
8	10		LH-729120	LH-747321	
	100		LH-729130	-	
	120		-	LH-747341	
	300		LH-729140	LH-747361	
	1,200		-	LH-747391	
12	10		LH-729220	LH-747421	
	100		LH-729230	-	
	120		-	LH-747441	
	300		LH-729240	LH-747461	
	1,200		-	LH-747491	

Event 2.  Tacta multi pack 할인	Product	Channel-Volume Range ( $\mu\ell$ )	Order no.
	Tacta 3-pack 10	1ch-10/100/1,000	LH-729670
	Tacta 3-pack 20	1ch-20/200/1,000	LH-729671
	Tacta 4-pack 20	1ch-10/20/200/1,000	LH-729672
	Tacta 4-pack 100	1ch-10/100/200/1,000	LH-729673
	Tacta 5-pack 10	1ch-10/100/200/1,000/5,000	LH-729674
	Tacta 5-pack 20	1ch-20/100/200/1,000/5,000	LH-729675
	Tacta 3+1-pack	1ch-10/200/1,000, 8ch-300	LH-729676
	Tacta 3-pack Max	1ch-1,000/5,000/10,000	LH-729677

QR  
CODE

Find out more