

Vivaspin Turbo RC 재 출시 기념 프로모션

QR
CODE

Vivaspin Turbo RC (48pc) 1박스 구매 시,
수동 피펫(Tacta) 1자루 무상 증정
기간 : 2023.12.31 까지



Features of Vivaspin Turbo RC

- 단시간에 Ultrafiltration & Diafiltration 가능
- 완벽한 Recovery와 최소 Dead Stop Volume
- 단단한 Housing으로 인한 최고의 내구성

프로모션 대상 제품

Order No.	Molecular Weight Cut Off
VS15TR12	5 kDa
VS15TR02	10 kDa
VS15TR22	30 kDa
VS15TR32	50 kDa
VS15TR42	100 kDa

※ 모든 제품에 대한 Sample Test 지원



수동 피펫(Tacta) 증정 품목



Volume Range(μl)

0.1 - 3

0.5 - 10

2 - 20

10 - 100

20 - 200

100 - 1000

500 - 5000

1 ml - 10ml

명함

싸토리우스코리아바이오텍(유)
경기도 성남시 분당구 판교역로 220 쏠리드스페이스 8층

E . marketing@sartorius.co.kr

www.sartorius.co.kr

무상증정(약 42만원 상당)

※위 품목 중 한 개 선택

Technical Specifications



	<i>Vivaspin 500</i>	<i>Vivaspin 2</i>	<i>Vivaspin Turbo 4</i>	<i>Vivaspin 6</i>	<i>Vivaspin 15R</i>	<i>Vivaspin Turbo15</i>	<i>Vivaspin 20</i>
Membrane area	0.5cm ²	1.2cm ²	3.2cm ²	2.5cm ²	3.9cm ²	7.2cm ²	6.0cm ²
Membrane	PES	PES,CTA, HY	PES	PES	HY	PES, RC	PES
Sample vol.	0.1 - 0.5ml	0.4 - 2ml	1 - 4ml	2 - 6ml	2 - 15ml	4 - 15ml	5 - 20ml

Ordering Information

MWCO	<i>Vivaspin 500</i>	<i>Vivaspin 2</i>	<i>Vivaspin Turbo 4</i>	<i>Vivaspin 6</i>	<i>Vivaspin 15R</i>	<i>Vivaspin Turbo15</i>	<i>Vivaspin 20</i>	
	PES	PES	PES	PES	HY	PES	RC	PES
2,000	-	-	-	-	VS15RH91	-	-	-
3,000	VS0191	VS0291	VS04T91	VS0691	-	VS15T91	-	VS2091
5,000	VS0111	VS0211	VS04T11	VS0611	VS15RH11	VS15T11	VS15TR11	VS2011
10,000	VS0101	VS0201	VS04T01	VS0601	VS15RH01	VS15T01	VS15TR01	VS2001
30,000	VS0121	VS0221	VS04T21	VS0621	VS15RH21	VS15T21	VS15TR21	VS2021
50,000	VS0131	VS0231	VS04T31	VS0631	-	VS15T31	VS15TR31	VS2031
100,000	VS0141	VS0241	VS04T41	VS0641	-	VS15T41	VS15TR41	VS2041
300,000	VS0151	VS0251	-	VS0651	-	-	-	VS2051
1,000,000	VS0161	VS0261	-	VS0661	-	-	-	VS2061
0.2µm	VS0171	VS0271	-	VS0671	-	-	-	VS2071



Find out more