

SARTORIUS

Minisart Best Seller Promotion

Minisart<Syringe Filter> 프로모션

기간 : 2023.10.31 까지



First in Class **Minisart<Syringe Filter>** 베스트셀러 TOP 10!

GET 25% OFF
Super SALE



※ 베스트셀러 TOP 10 정보확인 다음 장으로 ▶GO▶

QR
CODE

Find out more

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

Product Information

■ Minisart Top 10 LIST

Order No.	Membrane	Housing	Pore Size	Connector Outlet	Sterile	Ø in mm	Qty/Pk
S6555-FMOSK	SFCA	MBS	0.45µm	Male Luer Lock	Yes	28mm	50
S6534-FMOSK	SFCA	MBS	0.2µm	Male Luer Lock	Yes	28mm	50
17765-Q	RC	PP	0.45µm	Male Luer Slip	No	25mm	500
17821-K	RC	PP	0.2µm	Male Luer Slip	No	4mm	50
17762-Q	RC	PP	0.45µm	Male Luer Slip	No	15mm	500
17593-K	CA	MBS	1.2µm	Male Luer Lock	Yes	28mm	50
S7594-FMOSK	CA	MBS	5µm	Male Luer Lock	Yes	28mm	50
16532-K	PES	PP	0.22µm	Male Luer Lock	Yes	28mm	50
16541-K	PES	PP	0.22µm	Male Luer Slip	Yes	28mm	50
17575-K	PTFE	PP	0.2µm	Male Luer Slip	No	25mm	50

명함