

SARTORIUS



Vivaflow® SU

Ready, Set, Flow

Experience the unique simplicity and ultimate performance of Vivaflow® SU. Re-designed from the ground up, these next generation, laboratory-dedicated tangential flow filtration (TFF) cassettes offer:

- Unparalleled ease of use with intuitive setup and no optimization effort
- Economical running on standard pumps without the need for cassette holders
- Efficient concentration and buffer exchange for feed volumes from 0.1 - 1 L
- Reliable processing and high recoveries for more molecule types than ever before
- Maintained sample integrity without risk of carryover, in a safer working environment
- Increased sustainability with less single-use waste and more recyclable packaging

Ordering Information

Membrane Material	MWCO	Recommended Feed Volume	
		100 – 500 mL	500 – 1,000 mL
Hydrosart® regenerated cellulose (RC)	2 kDa	VF-S050H0002-IV	VF-S050H0002-SV
	5 kDa	VF-S050H0005-IV	VF-S050H0005-SV
	10 kDa	VF-S050H0010-IV	VF-S050H0010-SV
	30 kDa	VF-S050H0030-IV	VF-S050H0030-SV
	100 kDa	VF-S050H0100-IV	VF-S050H0100-SV
	300 kDa	VF-S050H0300-IV	VF-S050H0300-SV
Polyethersulfone (PES)	5 kDa	VF-S050P0005-IV	VF-S050P0005-SV
	10 kDa	VF-S050P0010-IV	VF-S050P0010-SV
	30 kDa	VF-S050P0030-IV	VF-S050P0030-SV
	50 kDa	VF-S050P0050-IV	VF-S050P0050-SV
	100 kDa	VF-S050P0100-IV	VF-S050P0100-SV
	300 kDa	VF-S050P0300-IV	VF-S050P0300-SV
	1,000 kDa	VF-S050P1000-IV	VF-S050P1000-SV
	0.2 µm	VF-S050P2000-IV	VF-S050P2000-SV

Equipment and Accessories

Description	Pack Size	Order No.
Peristaltic pump	1	VF-APD0001-1
Peristaltic pump head for 1.6 mm tubing	1	VF-APH0001-1
Cassette stand	1	VFA016
Diafiltration tubing	1	VF-ATD0001-1
Feed reservoir	1	VFA006

For More Information, Please Contact your Sales Representative