

SARTORIUS



Vivaflow[®]

Intuitive, Effortless
and Sustainable TFF
for the Lab

Groundbreaking Vivaflow[®] cassettes from Sartorius simplify TFF for researchers who value efficiency over complexity.

The ready-to-use cassettes are available in reusable or brand-new single use formats. In addition the modular design and comprehensive selection of membranes ensures you can always find the right choice to concentrate and buffer exchange any 0.1 – 5 L sample.



Plug and Play

Simple switch from centrifugal UF. Fast and intuitive set up right out of the box.



No Added Costs

Push your budget further. Built to run without expensive systems and holders.



Pre-Optimized Performance

Put down the calculator. Efficient UF | DF without process optimization.



Options for Any Sample

Ultimate Flexibility. Eliminate molecule loss with modular single-use or reusable cassettes.



Eco-Conscious Single Use

Reduce your research footprint. Vivaflow® SU uses 30.5 % less plastic and better packaging.

Ordering Information

Membrane Material	MWCO	Vivaflow® SU ¹	Vivaflow® SU ¹	Vivaflow® 50R ²	Vivaflow® 200 ²
Hydrosart® regenerated cellulose (RC)	2 kDa	VF-S050H0002-IV	VF-S050H0002-SV	-	VF20H9
	5 kDa	VF-S050H0005-IV	VF-S050H0005-SV	VF05H1	VF20H1
	10 kDa	VF-S050H0010-IV	VF-S050H0010-SV	VF05H0	VF20H0
	30 kDa	VF-S050H0030-IV	VF-S050H0030-SV	VF05H2	VF20H2
	100 kDa	VF-S050H0100-IV	VF-S050H0100-SV	VF05H4	VS20H4
	300 kDa	VF-S050H0300-IV	VF-S050H0300-SV	-	-
Polyethersulfone (PES)	3 kDa	-	-	-	VF20P9
	5 kDa	VF-S050P0005-IV	VF-S050P0005-SV	-	VF20P1
	10 kDa	VF-S050P0010-IV	VF-S050P0010-SV	-	VF20P0
	30 kDa	VF-S050P0030-IV	VF-S050P0030-SV	-	VF20P2
	50 kDa	VF-S050P0050-IV	VF-S050P0050-SV	-	VF20P3
	100 kDa	VF-S050P0100-IV	VF-S050P0100-SV	-	VF20P4
	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	-	VF20P7	

¹ Order numbers ending -IV include two cassettes and two tubing kits to run each cassette individually. Order numbers ending -SV include two cassettes and one tubing kit to run both cassettes in series.

² Includes one cassette and one tubing kit. To run two cassettes in parallel, additional accessories are required and can be ordered separately.

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
Peristaltic pump head for 2.4 mm tubing	1	VF-APH0002-1
Cassette stand	1	VF-AST0001-1
Feed reservoir	1	VF-ARV0500-1
Flow restrictors, 0.4 and 0.8 mm	2	VF-AC-0002-V
Tubing for diafiltration	1	VF-ATD0001-1
Tubing for individual cassettes	1	VF-ATI0011-1
Tubing for 2x 50 cm ² cassettes in parallel	1	VF-ATP0011-1
Tubing for 2x 200 cm ² cassettes in parallel	1	VF-ATP0012-1

For More Information, Please Contact your Sales Representative