

Vivapore®

Keep Still and Concentrate



Product Information

Vivapore® solvent absorption ultrafilters for general laboratory use offer a reliable and user-friendly alternative for sample concentration. They are particularly suitable for concentrating proteins with molecular weights ≥ 25 kDa in research and clinical laboratories.

Although centrifugal ultrafilters have become the gold standard for rapid concentration of macromolecules from small process volumes, they are not always the most practical choice.

In cases where access to centrifuges is limited or their use is impractical (e.g. in laboratories processing high quantities of samples on a daily basis), Vivapore® enables sample concentration without the need for additional equipment.

Equipped with high-performance polyethersulfone (PES) membranes, Vivapore® units also feature a compact footprint and at-a-glance process monitoring, while eliminating the risk of sample loss and cross-contamination.

Features

Concentrate Equipment-Free

Solvent absorption is ideal where centrifuges are not available and for high throughput applications.

Fill and Go

Simply add your sample and leave it to reach the desired final volume, with no risk of concentration to dryness.

Tailored Process Monitoring

Printed markings help to reduce the guesswork when each sample needs to reach its own specific concentration factor.

Maintained Sample Integrity

Self-contained, single use units ensure contamination-free concentration of individual samples.

Sensitive Molecule Friendly

Benchtop or refrigerated low pressure processing for shear and temperature sensitive molecules.

Expand or Exchange¹

Optional reservoirs provide the flexibility to double sample capacity or concentrate and desalt in one step.

¹ Expansion reservoirs are available for Vivapore® 10 only.

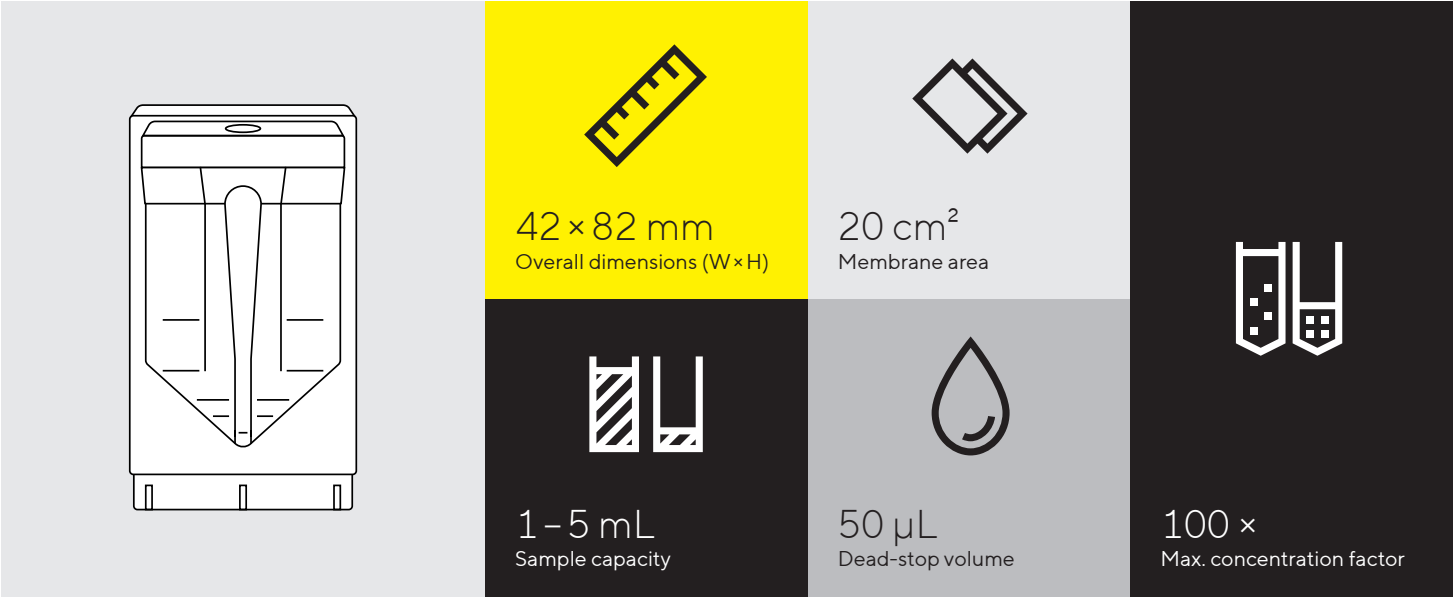
Applications

- Concentrating samples prior to *in vitro* diagnostics
- Concentrating shear-sensitive molecules
- Concentrating temperature-sensitive samples
- Desalting to reduce interference in downstream analysis

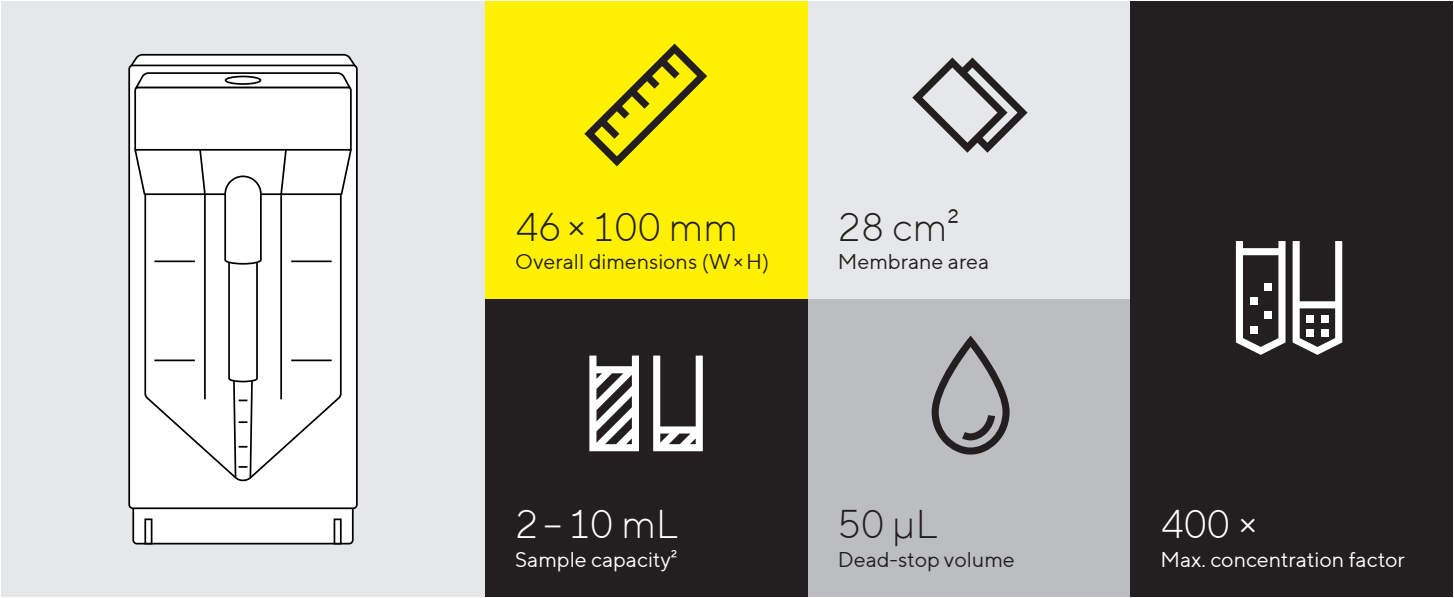


Technical Specifications

Vivapore® 5



Vivapore® 10



² Up to 20 mL sample capacity when used with an expansion reservoir.

Materials

Housing	Styrene acrylonitrile (SAN)
Membrane	Polyethersulfone (PES)
Absorbent	Cellulose
Stand	Acrylonitrile butadiene styrene (ABS)
Reservoir	Polypropylene (PP)
Packaging	Cardboard (PAP) Polyethylene (LDPE)

Typical Performance

Typical process time and protein recovery for 5, 10 and 20 mL starting volumes (Vivapore® 5, Vivapore® 10 and Vivapore® 10 with expansion reservoir, respectively) at 20 °C.

Product	Protein	MW	10× Concentration Factor		50× Concentration Factor	
			Time	Recovery	Time	Recovery
Vivapore® 5	Cytochrome c	12.6 kDa	35 min	90%	65 min	91%
	BSA	66 kDa	30 min	92%	45 min	90%
	IgG	150 kDa	40 min	75%	50 min	53%
Vivapore® 10	Cytochrome c	12.6 kDa	75 min	90%	70 min	88%
	BSA	66 kDa	55 min	92%	50 min	90%
	IgG	150 kDa	70 min	77%	65 min	65%
Vivapore® 10 with reservoir	Cytochrome c	12.6 kDa	150 min	92%	160 min	90%
	BSA	66 kDa	115 min	92%	105 min	92%
	IgG	150 kDa	160 min	78%	140 min	74%

Ordering Information

Description	Package Contents	Order No.
Vivapore® 5 7.5 kDa MWCO PES	30 units	VP-S005P0008-3
	1 stand	
	1 quick start guide	
Vivapore® 10 7.5 kDa MWCO PES	30 units	VP-S010P0008-3
	1 stand	
	1 expansion reservoir	
	1 quick start guide	
Stand	6 units	VP-AST0001-C
Expansion reservoir ³	10 units	VP-ARV0010-D

³ Expansion reservoirs are compatible with Vivapore® 10 only.

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further information, visit
www.sartorius.com