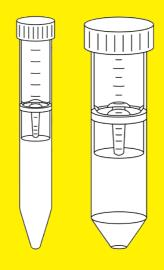
Instructions for Use

Vivaspin® 6 | 20

Centrifugal Ultrafiltration Units for General Laboratory Use



3104655-001-00





Contents

1	Abo	ut thes	e Instructions	5
	1.1	Scope	9	5
	1.2	Targe	t Groups	6
	1.3	Symb	ols Used	7
		1.3.1	Warnings in Operation Descriptions	7
		1.3.2	·	
2	Safe	ty Instr	ructions	8
	2.1	Gene	ral functions	8
	2.2	Perso	nnel Qualification	9
	2.3	Signif	icance of these Instructions	9
	2.4	Funct	ionality of the Product	9
3	Proc	luct De	escription	10
	3.1	Vivası	oin [®] 6	10
	3.2	Vivası	pin [®] 20	11
	3.3	Produ	uct Symbols	12
4	Proc	ess Pre	eparation	13
	4.1	Scope	e of Delivery	13
	4.2	Unpa	cking	13
5	Оре	ration.		14
	5.1	Filtrat	ion in Centrifuges	14
		5.1.1	Pre-Rinsing the Product	14
		5.1.2	Sanitizing the Product	14
		5.1.3	Applying the Sample	15
		5.1.4	Inserting the Product into the Centrifuge	16
		5.1.5	Performing Filtration	16
		5.1.6	Removing the Sample	16
		5.1.7	Removing the Concentrator from the Filtrate Container.	17
	5.2	Filtrat	ion using Gas Pressure	18
	5.3	Conti	nuous Diafiltration (Vivaspin® 20 only)	19
	5.4	Desal	ting or Buffer Exchange	20

Contents

6	Stor	age		21
	6.1	Storin	ng the Product	21
7	Disp	osal		22
	7.1		ntaminating the Product	
	7.2		osing of the Product	
8	Tech	nical S	pecifications	23
	8.1		nsions	
	8.2		rials	
	8.3		ent Conditions	
	8.4		ating conditions	
		8.4.1	Filtration Volumes	
		8.4.2	Centrifugation Limit Values	
			Pressurization Limit Values	
	8.5		oment Required	
		8.5.1	Pipettes	
		8.5.2		
		8.5.3	Using gas pressure	
	8.6		zing Methods	
	8.7		nical Compatibility	
	8.8		al Performance Characteristics	
	0.0	8.8.1	Vivaspin [®] 6	
			Vivaspin® 20	
			r	
9	Orde	ering In	nformation	31

1 About these Instructions

1.1 Scope

These instructions are part of the product. These instructions apply to the following versions of the product:

Vivaspin® 6	Quantity	Prod. no.
3 kDa	25 100	VS0691 VS0692
5 kDa	25 100	VS0611 VS0612
10 kDa	25 100	VS0601 VS0602
30 kDa	25 100	VS0621 VS0622
50 kDa	25 100	VS0631 VS0632
100 kDa	25 100	VS0641 VS0642
300 kDa	25 100	VS0651 VS0652
1,000 kDa	25 100	VS0661 VS0662
0.2 μm	25 100	VS0671 VS0672

Vivaspin® 20	Quantity	Prod. no.
3 kDa	12 48	VS2091 VS2092
5 kDa	12 48	VS2011 VS2012
10 kDa	12 48	VS2001 VS2002
30 kDa	12 48	VS2021 VS2022
50 kDa	12 48	VS2031 VS2032
100 kDa	12 48	VS2041 VS2042
300 kDa	12 48	VS2051 VS2052
1,000 kDa	12 48	VS2061 VS2062
0.2 μm	12 48	VS2071 VS2072

1.2 Target Groups

The instructions are designed for the following target groups. The target groups must possess the knowledge listed below.

Target Group	Knowledge and Qualifications		
User	The user is familiar with how to use the product and with the associated work processes. The user is familiar with the hazards that may arise when working with the product, and is able to avoid these hazards. The user has received training on how to use the product.		
Operator	The operator of the product is responsible for compliance with safety requirements and workplace safety regulations. The operator must ensure that anyone working with the product has access to the relevant information and is trained to work with the product.		

1.3 Symbols Used

1.3.1 Warnings in Operation Descriptions

NOTICE

Denotes a hazard that may result in property damage if it is **not** avoided.

1.3.2 Other Symbols

- Required action: Describes actions that must be carried out.

 The actions in the sequence must be carried out in succession.
- Result: Describes the result of the actions carried out.

2 Safety Instructions

2.1 General functions

The product is intended for the ultrafiltration and | or diafiltration of biological and aqueous solutions. The sample solutions and volumes used must be suitable for the product.

The filtration process must be carried out in a centrifuge. Macromolecules that are sufficiently larger than the nominal pore size of the membrane are retained above the membrane and progressively concentrated. The vertical membrane inhibits membrane fouling while the built-in dead stop impedes concentration to dryness and loss of sample.

For Vivaspin® 20, the filtration process may alternatively be carried out by using gas pressure as well as by using a diafiltration cup. It is also possible to combine gas pressure with centrifugation (pressure-fugation).

If Vivaspin® 20 is used with a diafiltration cup: During filtration salts and other contaminating micromolecules can be removed from the sample by continous diafiltration.

The product is supplied non-sterile. It is intended for single use and must be disposed of after one use.

The product is intended exclusively for use in accordance with these instructions. Any further use beyond this is considered improper.

Operating Conditions for the Product

The product is intended for general laboratory use.

The product may only be used with the equipment and under the operating conditions described in the Technical Data section of these instructions.

2.2 Personnel Qualification

Persons without sufficient knowledge in the safe use of the device can injure themselves and others.

If a specific qualification is required for an activity: The target group is indicated. If no qualification is specified: The activity can be carried out by the target group "Operator".

2.3 Significance of these Instructions

Failure to follow the instructions might have serious consequences, e.g. danger to individuals.

- ▶ Read the instructions carefully and completely. The instructions for action build on each other.
- ► Ensure that the information contained in these instructions is available to all individuals working with the product.

2.4 Functionality of the Product

A damaged product or worn parts can lead to malfunctions or cause hazards which are difficult to identify.

▶ Only operate the product when it is safe and in perfect working order.

3 Product Description

3.1 Vivaspin® 6

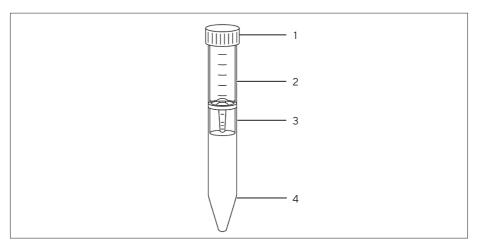


Fig. 1: Product overview (example)

Pos.	Description
1	Concentrator cap
2	Concentrator
3	Twin vertical membranes
4	Filtrate container

3.2 Vivaspin® 20

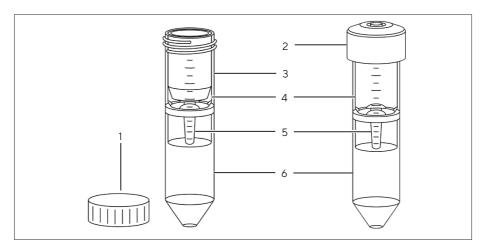


Fig. 2: Product overview (example)

Pos.	Description
1	Concentrator cap
2	Pressure cap (optional accessory, VCA200)
3	Diafiltration cup (optional accessory, VSA005)
4	Concentrator
5	Twin vertical membranes
6	Filtrate container

3.3 Product Symbols

Symbol	Definition
REF	Catalogue number
②	Do not reuse
\square	Use by
LOT	Batch code
	Manufacturer
X	Temperature limitation
MON	Non-sterile product

4 Process Preparation

4.1 Scope of Delivery

Article	Qty	
Product packed in a cardboard box		
Vivaspin® 6	25 or 100	
Vivaspin® 20	12 or 48	
Instructions for Use	1	

4.2 Unpacking

- ▶ **NOTICE** Risk of product malfunctions due to exceeding the usability! Check the usability of the product (see specification on packaging). Dispose of products for which the usability has been exceeded.
- ► Unpack the product.

5 Operation

5.1 Filtration in Centrifuges

5.1.1 Pre-Rinsing the Product

Membranes in the product may contain traces of glycerin. If this substance can interfere with the analysis of the sample: The membranes may be rinsed before filtration.

Procedure

- ▶ Remove the concentrator cap.
- ▶ Use a pipette to apply a filling volume of buffer solution or deionized water into the concentrator.
- ▶ Replace the concentrator cap.
- ► Wash the buffer solution or deionized water through the membranes by centrifugation.
- ▶ Empty the concentrator and filtrate container.
- ▶ If the pre-rinsed product is not used immediately: Cover the surface of the membrane with buffer solution or water and store the product in the refrigerator. The membranes must not dry out.

5.1.2 Sanitizing the Product

The product can be sanitized before use. The sanitizing method must be suitable for the product (see Chapter "8.6 Sanitizing Methods", page 26).

- Remove the concentrator cap.
- ▶ Sanitize the product using the desired sanitizing method.
- Empty the product.

5.1.3 Applying the Sample

It is recommended that a pipette is used to apply the sample into the product. The pipette must be compatible with the product (see Chapter "8.5.1 Pipettes", page 26).

Please ensure that the molecular weight cut-off (MWCO) of the product is suitable for the size of the target molecule to be concentrated. In order to ensure maximum recovery of the target molecule, it is recommended to select a MWCO that is at least 50% below the size of the target molecule.

NOTICE

Risk of product malfunctions due to using unsuitable samples!

▶ Only pour suitable samples into the product (see Chapter "8.7 Chemical Compatibility", page 27).

NOTICE

Risk of product malfunctions or damage to the centrifuge due to exceeding the maximum filling volume!

▶ Do **not** exceed the maximum filling volume (see Chapter "8.4 Operating conditions", page 24).

- ▶ Check whether the MWCO of the product is suitable for the application.
- ► Remove the concentrator cap.
- ▶ Apply the sample into the product using a pipette. Comply with the maximum filling volume.
- Replace the concentrator cap.

5.1.4 Inserting the Product into the Centrifuge

Procedure

- ▶ NOTICE Risk of product malfunctions or damage to the centrifuge! Only use the product in suitable centrifuges (see Chapter "8.5.2 Centrifuges", page 26).
- Insert the product into the centrifuge.
- ▶ If a centrifuge with fixed-angle rotor is used: Place the product into the centrifuge so that the printed volume graduations on the concentrator are facing upwards.

5.1.5 Performing Filtration

Procedure

- ▶ **NOTICE** Risk of product malfunctions or damage to the centrifuge. Comply with the approved centrifugation limit values (see Chapter "8.4.2 Centrifugation Limit Values", page 25).
- ► Centrifuge the product in the centrifuge until the desired concentration level is achieved.

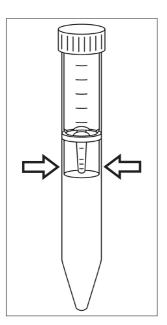
5.1.6 Removing the Sample

- ▶ If the filtration or concentration is complete: Remove the product from the centrifuge.
- ► Remove the concentrator cap.
- Recover the sample from dead stop pocket of the concentrator using a pipette.
- ▶ If the membrane was pre-rinsed before filtration: Decant the filtrate and concentrate.

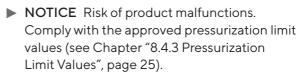
5.1.7 Removing the Concentrator from the Filtrate Container

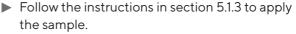


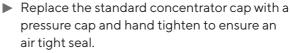
➤ To release the tube from the concentrator, pinch the tube on either side of the printed volume graduations – to press it into an oval shape – before removing it with a twisting action.



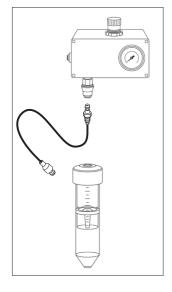
5.2 Filtration using Gas Pressure



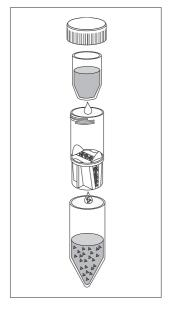




- When using the Air Pressure Controller (APC): Replace the female coupling on the extension line with the appropriate charge valve (VCAO05) and insert this into the inlet valve of the pressure cap to pressurize the product.
- ➤ Stand the product in a centrifuge tube rack until the sample is concentrated to the desired level. Alternatively, insert the product into a centrifuge and centrifuge the product until the desired concentration level is achieved.
- ▶ If the filtration or concentration is complete: De-pressurize the product by unscrewing the pressure cap.
- ► Recover the sample from dead stop pocket of the concentrator using a pipette.



5.3 Continuous Diafiltration (Vivaspin® 20 only)



- Apply up to 2 mL of sample into the concentrator. If the sample for diafiltration has a larger volume: Concentrate the sample to 2 mL and decant the filtrate.
- ► Insert a diafiltration cup into the concentrator and fill it with 10 mL buffer solution.
- ► Replace the concentrator cap.
- Insert the product into the centrifuge and centrifuge the product until diafiltration is complete.
- Salts and contaminants are progressively cleared through membrane and into filtrate vessel.
- ▶ Remove the product from the centrifuge.
- ► Remove the concentrator cap and the diafiltration cup.
- Recover the sample from dead stop pocket of the concentrator using a pipette.

5.4 Desalting or Buffer Exchange

- ► Concentrate the sample to the desired level.
- ▶ Remove the concentrator cap.
- Discard the filtrate.
- ► Refill the concentrator with an appropriate exchange buffer.
- ► Concentrate the sample again.
- ➤ Repeat the process until the original buffer and | or contaminating microsolute has been sufficiently removed.
- ► If the desalting or buffer exchange is complete: Recover the sample.

6 Storage

6.1 Storing the Product

If the product has been unpacked and membranes have been pre-rinsed: The membranes must be protected against drying out. For this purpose, the membranes must be stored in a moist and cool condition.

NOTICE

Risk of damage to the product due to improper storage!

Comply with the storage specifications.

- If the product is packaged: Store the product in the packaging.
- ▶ If the product has been unpacked and the membranes have been prerinsed:
 - ▶ Remove the concentrator cap.
 - ► Cover the membranes with buffer solution or water.
 - ► Replace the concentrator cap.
- ➤ Store the product according to the ambient conditions (see Chapter "8.3 Ambient Conditions", page 24).

7 Disposal

7.1 Decontaminating the Product

If the product has come into contact with hazardous substances: Steps must be taken to ensure proper decontamination and declaration. The operator of the product is responsible for adhering to local government regulations on the proper decontamination and declaration for transport and disposal.

Procedure

► If the product has come into contact with hazardous substances: Decontaminate the product.

7.2 Disposing of the Product

The product must be disposed of properly. The packaging is made of environmentally friendly materials that can be used as secondary raw materials.

Requirements

The product must be decontaminated.

- ▶ Dispose of the product in accordance with local government regulations.
- ▶ Dispose of the packaging in accordance with local government regulations.

8 Technical Specifications

8.1 Dimensions

	Vivaspin® 6		Vivaspin® 20	
	Unit	Value	Unit	Value
Length x Diameter	mm	122 x 17	mm	116 x 30 125 x 30 (with pressure cap)
Active membrane surface	cm²	2.5	cm²	6.0
Weight	g	10	g	25

8.2 Materials

	Vivaspin® 6	Vivaspin® 20
Concentrator Filtrate container	Polycarbonate	Polycarbonate
Concentrator cap	Polypropylene	Polypropylene
Pressure cap	-	Polyoxymethylene/ Aluminium
Membrane	Polyethersulfone	Polyethersulfone

8.3 Ambient Conditions

	Unit	Value
Storage temperature		
When packed	°C	+15 - +30
When unpacked, with membranes kept moist	°C	+2 - +8

8.4 Operating conditions

8.4.1 Filtration Volumes

		Centrifuge with swing bucket rotor	Centrifuge with fixed-angle rotor (25°)
	Unit	Value	Value
Vivaspin® 6			
Filling volume, maximum	mL	6	6
Membrane hold-up volume, minimum	μL	< 10	< 10
Dead stop volume ¹	μL	30	30
Vivaspin® 20			
Filling volume, maximum	mL	20	14
Filling volume with pressure cap, maximum	mL	15	-
Membrane hold-up volume, minimum	μL	< 20	< 20
Dead stop volume ¹	μL	50	50

 $^{^1\}text{The}$ dead stop volume may vary depending on the type and concentration of the sample, operating temperature and | or centrifuge rotor

8.4.2 Centrifugation Limit Values

		Centrifuge with swing bucket rotor	Centrifuge with fixed angle rotor
	Unit	Value	Value
Vivaspin® 6			
3 - 50 kDa PES	g	4,000	8,000
100 kDa - 0.2 μm PES	g	4,000	6,000
Vivaspin® 20			
3 - 50 kDa PES	g	4,000	6,000
100 kDa - 0.2 μm PES	g	3,000	6,000

8.4.3 Pressurization Limit Values

		Pressure	Pressure-Fuge with swing bucket rotor
	Unit	Value	Value
Vivaspin® 20			
3 - 50 kDa PES	bar	5	5
	g	-	3,000
100 kDa- 0.2 µm PES	bar	5	5
	g	-	2,000

8.5 Equipment Required

8.5.1 Pipettes

Pasteur pipette, variable volume or fixed volume pipette for sample application and concentrate or filtrate retrieval.

8.5.2 Centrifuges

Centrifuge with swing bucket rotor or fixed-angle rotor (minimum 25°) accepting conical base tubes.

	Required	Required carriers		
	Unit	Value		
Vivaspin® 6				
Volume	mL	15		
Diameter	mm	17		
Vivaspin® 20				
Volume	mL	50		
Diameter	mm	30		

8.5.3 Using gas pressure

	Vivaspin [®] 20
Approved	- Pressure cap
devices	 Charge valve for pressure cap
	 Air Pressure Controller or suitable pressure regulator

8.6 Sanitizing Methods

Rinsing with 70% ethanol or with sanitizing gas mixture, e.g. ethylene oxide

Not suitable for autoclaving

8.7 Chemical Compatibility

Chemical, biological and aqueous solutions with appropriate compatibility for the materials of the product (2 hr contact time)

Examples of compatible chemical solutions				
Acetic Acid (25%)	Lactic Acid (5%)			
Acetone (10%)	Mercaptoethanol (10 mM)			
Acetonitrile (10%)	Methanol (60%)			
Ammonium Hydroxide (5%)	Nitric Acid (10%)			
Ammonium Sulphate (saturated)	Phenol (1%)			
Benzene (100%)	Phosphate Buffer (1 M)			
n-Butanol (70%)	Polyethylene Glycol (10%)			
Chloroform (1%)	Pyridine (100%)			
Dimethyl Formamide (10%)	Sodium Carbonate (20%)			
Dimethyl Sulfoxide (5%)	Sodium deoxychloate (5%)			
Ethanol (70%)	Sodium dodecylsulfate (0.1 M)			
Ethyl Acetate (100%)	Sodium Hydroxide (2 M)			
Formaldehyde (30%)	Sodium Hypochlorite (200 ppm)			
Formic acid (5%)	Sodium Nitrate (1%)			
Glycerine (70%)	Sulfamic Acid (5%)			
Guanidine HCI (6 M)	Tetrahydrofuran (5%)			
Hydrocarbons, aromatic	Trifluoroacetic Acid (10%)			
Hydrocarbons, chlorinated	Triton®* X-100 (0.1%)			
Hydrochloric Acid (1 M)	Tween®** 20 (0.1%)			
Imidazole (500 mM)	Urea (8 M)			
Isopropanol (70%)				

^{*} Triton® is a registered trademark of Union Carbide Corp.

^{**} Tween® is a registered trademark of ICI Americas Inc.

8.8 Typical Performance Characteristics

8.8.1 Vivaspin® 6

Swing bucke	·	to 30x at 20°			
6 mL		i inca arigie	Fixed angle (25°)		
		6 mL			
Time (min)	Solute Recovery	Time (min)	Solute Recovery		
nL (12.4 kDa)	1				
-	-	90	97%		
20	98%	12	98%		
13	98%	10	98%		
12	98%	9	97%		
a)					
18	96%	15	95%		
17	96%	14	95%		
15	91%	12	91%		
MEM + 10%	FCS (0.055 _k	ım)			
-	-	25	99%		
Latex beads 0.004% in DMEM + 10% FCS (0.24 μm)					
-	-	4	99%		
visiae)					
4	97%	3	97%		
	nL (12.4 kDa) - 20 13 12 a) 18 17 15 MEM + 10% - WEM + 10% - visiae)	Recovery nL (12.4 kDa) 20 98% 13 98% 12 98% a) 18 96% 17 96% 15 91% MEM + 10% FCS (0.055 µ MEM + 10% FCS (0.24 µn visiae)	Recovery nL (12.4 kDa) 90 20 98% 12 13 98% 10 12 98% 9 a) 18 96% 15 17 96% 14 15 91% 12 MEM+10% FCS (0.055 µm) - 25 MEM+10% FCS (0.24 µm) - 4 visiae)		

8.8.2 Vivaspin® 20

Time to concentrate up to 30x at 20°C								
 Mode				<u>'</u>	Bench top		Pressure-fuge	
Rotor		Centrifuge Centrifuge Swing bucket Fixed angle (25°)			Pressure		Swing bucket	
Start volume	20 mL		14 mL		10 mL		10 mL	
	Time (min)	Solute Re- covery	Time (min)	Solute Re- covery	Time (min)	Solute Re- covery	Time (min)	Solute Re- covery
Cytochrome	c 0.25 m	ng/mL (1	2:4 kDa))				
3 kDa PES	110	97%	180	96%	60	96%	-	-
BSA 1.0 mg/m	nL (66 k	Da)						
5 kDa PES	23	99%	29	99%	50	98%	14	98%
10 kDa PES	16	98%	17	98%	32	97%	8	97%
30 kDa PES	13	98%	15	98%	32	97%	8	97%
IgG 0.25 mg/	mL (160	kDa)						
30 kDa PES	27	97%	20	95%	46	94%	13	97%
50 kDa PES	27	96%	22	95%	46	93%	13	96%
100 kDa PES	25	91%	20	90%	42	88%	12	94%

	Time to	concer	ntrate u _l	o to 30x	at 20°C	;			
Mode	Centrif	fuge Centrifuge		uge	Bench top		Pressure-fuge		
Rotor	Swing I	oucket	ucket Fixed an (25°)		Pressui	Pressure		Swing bucket	
Start volume	20 mL		14 mL		10 mL		10 mL		
	Time (min)	Solute Re- covery	Time (min)	Solute Re- covery	(min)	Solute Re- covery	(min)	Solute Re- covery	
Latex beads C	0.004%	in DMEI	M + 10%	FCS (0.	055 µm)			
30 kDa PES	20	99%	35	99%	10	99%	-	-	
Latex beads 0	0.004%	in DMEI	M + 10%	FCS (0.	24 µm)				
1,000 kDa PES	4	99%	12	99%	4	99%	-	-	
Yeast 1.0 mg/	mL (S. c	erevisia	e)						
0.2 µm PES	15	95%	5	95%	20	95%	2	95%	

9 Ordering Information

Vivaspin® 20 accessories	Quantity	Prod. no.
Air pressure controller (APC)	1	VCA002
Charge valve for pressure cap VCA200	1	VCA005
Diafiltration cups	12	VSA005
Female coupling	1	VCA010
Male coupling	1	VCA011
Replacement extension line	1	VCA012
Vivaspin® 20 pressure cap	1	VCA200

Sartorius Stedim Lab Ltd. Sperry Way, Stonehouse GL10 3UT, UK

Phone: +44 1453 821972 www.sartorius.com

The information and figures contained in these instructions correspond to the version date specified below.

Sartorius reserves the right to make changes to the technology, features, specifications and design of the equipment without notice.

Masculine or feminine forms are used to facilitate legibility in these instructions and always simultaneously denote all genders.

Copyright notice:

These instructions, including all components, are protected by copyright.

Any use beyond the limits of the copyright law is not permitted without our approval.

This applies in particular to reprinting, translation and editing irrespective of the type of media used.

Last updated: 02 | 2023

© 2023 Sartorius Stedim Lab Ltd. Sperry Way, Stonehouse GL10 3UT, UK KS | Publication No.: SLU6092-e230214 Material No.: 85030-511-57