



# PD MiniTrap™ G-25

## Instructions for Use

# Table of Contents

1	Introduction .....	3
2	Description .....	4
3	Advice on handling .....	6
4	Column assembly .....	7
5	Gravity protocol .....	8
6	Spin protocol .....	11
7	Recovery and desalting capacity .....	13
8	Column characteristics .....	14
9	Ordering information .....	15

# 1 Introduction

## Important

Read these instructions carefully before using the products.

## Intended use

The products are intended for research use only, and shall not be used in any clinical or *in vitro* procedures for diagnostic purposes.

## Safety

For use and handling of the products in a safe way, refer to the Safety Data Sheets.



### **CAUTION**

The column storage solution, 0.15% Kathon CG/ICP biocide, is potentially allergenic. Use gloves when discarding the storage solution.

## 2 Description

PD MiniTrap™ G-25 columns contain *Sephadex™ G-25 resin*, which allows rapid *group separation* of high molecular weight substances from low molecular weight substances.

PD MiniTrap G-25 columns are used for desalting, buffer exchange and sample clean up. Small molecules like salt, free labels and other impurities are efficiently separated from the high molecular weight substances of interest. The chromatography technique is gel filtration and molecules are separated on the basis of differences in size.

- Molecules larger than the largest pores in the Sephadex matrix are excluded from the matrix and are eluted first, in or just after the void volume. The void volume is the column volume outside the Sephadex matrix.
- Molecules smaller than the largest pores in the Sephadex matrix will penetrate the pores to varying extent. They have a larger accessible column volume than the large molecules and therefore they elute after the large molecules just before one total column volume of buffer has passed through the column.

Group separation can be made using two different protocols, gravity protocol and spin protocol.

Cytiva provides an assortment of sample clean-up products. The different formats available are summarized in [Table 1, on page 5](#).

**Table 1.** Product overview

Clean Up product	Exclusion limit, M <sub>r</sub>	Bed volume	Sample volume gravity protocol <sup>1</sup>	Sample volume spin protocol <sup>1</sup>
PD SpinTrap™ G-25	5000	0.5 mL	-	100 to 180 µL
PD MultiTrap™ G-25	5000	0.5 mL	-	70 to 130 µL
PD MiniTrap G-25	5000	2.1 mL	0.1 to 0.5 mL	0.2 to 0.5 mL
PD MidiTrap™ G-25	5000	3.5 mL	0.5 to 1.0 mL	0.75 to 1.0 mL
PD-10 Desalting Columns	5000	8.3 mL	1.0 to 2.5 mL	1.75 to 2.5 mL
PD MiniTrap G-10	700	2.1 mL	0.1 to 0.3 mL	-
PD MidiTrap G-10	700	5.3 mL	0.4 to 1.0 mL	-

<sup>1</sup> Recommended sample volumes.

## 3 Advice on handling

### Protocol selection

The separation can be made using two different protocols, gravity protocol or spin protocol, see Table 2 below for protocol overview.

### Gravity protocol

The liquid passes through the column by gravity force.

- There is a slightly higher recovery and desalting capacity using gravity protocol compared to when using the spin protocol.
- The applied sample is diluted.

### Spin protocol

Additional gravity force is added by spinning the column in a centrifuge for some protocol steps.

- There is no dilution of the sample.

**Table 2.** Protocol overview

Protocol	Sample volume	Elution buffer	Dilution factor	Desalting capacity
Gravity	0.1 to 0.5 mL	1.0 mL	2 times <sup>1</sup>	>98%
Spin	0.2 to 0.5 mL	None	None	>90%

<sup>1</sup> Two times dilution valid if 0.5 mL sample volume is used.

### Recovery

The recovery of applied amount sample is dependent on type of protein or other biomolecule. Typically the recovery is in the range 70–90%. An increase in sample concentration can improve recovery.

## Equilibration

- It is critical to equilibrate the column since UV absorbing stabilizers are used in column packing.
- Equilibration is most conveniently made by gravity also when using the spin protocol.

## Sample Application

- The MiniTrap column is intended for sample volumes up to 0.5 mL.
- For sample volumes less than 0.5 mL, allow the sample to enter the packed bed completely and then add equilibration buffer (stacker volume) so that the total volume of sample and buffer added equals 0.5 mL.
- Allow the sample to enter the packed bed completely before any addition of buffer for elution.

## Elution

- For higher recoveries, use stacker volumes (spin protocol).

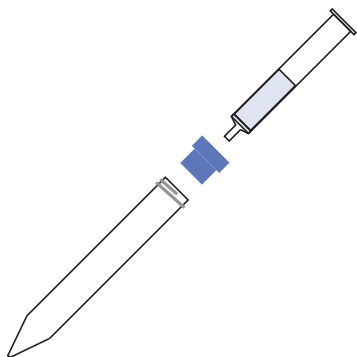
## Centrifugation

For better result if using a fixed angle rotor centrifuge: Put the columns in same direction in all centrifugation steps (spin protocol).

# 4 Column assembly

For use in the spin protocol the PD MiniTrap G-25 column must be assembled with adapter and collection tube as shown below in Figure 1.

**Note:** This column assembly may also be used for convenient handling of columns when using the gravity protocol.



**Fig 1.** Assembly of column, adapter and collection tube.

## 5 Gravity protocol

### PD MiniTrap G-25 preparation

Step	Action
------	--------

- |   |  |
|---|--|
| 1 | Remove the top cap and pour off the column storage solution. |
| 2 | Remove the bottom cap.                                       |



## Column equilibration

Step	Action
------	--------

- |   |   |
|---|---|
| 1 | Fill up the column with equilibration buffer and allow the equilibration buffer to enter the packed bed completely. |
| 2 | Repeat twice.   |
| 3 | Discard the flowthrough.  |

**Note:**

*In total, approximately 8 mL equilibration buffer must be used for all three steps.*

## Sample application

Step	Action
------	--------

- |   |   |
|---|---|
| 1 | Add maximum 0.5 mL of sample to the column.   |
| 2 | For sample volumes less than 0.5 mL, add equilibration buffer to adjust the volume up to 0.5 mL after the sample has entered the packed bed completely. |
| 3 | Let the sample or equilibration buffer enter the packed bed completely.   |
| 4 | Discard the flowthrough.  |

## Elution

Step	Action
------	--------

- |   |   |
|---|---|
| 1 | Place a test tube for sample collection under the column. |
|---|---|

<b>Step</b>	<b>Action</b>
-------------	---------------

---

- |          |  |
|----------|--|
| <b>2</b> | Elute with 1.0 mL buffer and collect the eluate.<br>A typical elution profile is shown in <a href="#">Fig. 2, on page 13</a> . |
|----------|--|
-

## 6 Spin protocol

### PD MiniTrap G-25 preparation

Step	Action
1	Remove the top cap and pour off the column storage solution.
2	Remove the top filter using forceps.
3	Remove the bottom cap.
4	Put the PD MiniTrap G-25 into a 15 mL collection tube by using the column adapter, see <a href="#">Fig. 1, on page 8</a> .

### Column equilibration

Step	Action
1	Fill up the column with equilibration buffer and allow the equilibration buffer to enter the packed bed completely.
2	Repeat once.
3	Discard the flow-through.
4	Fill up the column a third time with equilibration buffer and spin down at $1000 \times g$ for 2 minutes.
5	Discard the flow-through.

**Note:**  
*About 8 mL equilibration buffer should be used in total for all three steps.*

## Sample application

Step	Action
------	--------

---

- |   |  |
|---|--|
| 1 | Add sample (0.2 to 0.5 mL) slowly in the middle of the packed bed. |
|---|--|
- 

## Elution

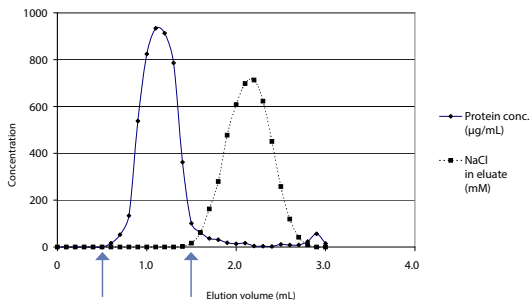
Step	Action
------	--------

---

- |   |  |
|---|--|
| 1 | Place the PD MiniTrap G-25 into a new 15 mL collection tube. |
| 2 | Elute by centrifugation $1000 \times g$ for 2 minutes.       |
| 3 | Collect the eluate.  |
-

## 7 Recovery and desalting capacity

The following experiment is included as an example of a desalting experiment using the gravity protocol. A PD MiniTrap G-25 column was equilibrated with Milli-Q water. 500  $\mu\text{L}$  of bovine serum albumin (1000  $\mu\text{g}/\text{mL}$ ) in 1 M NaCl was applied onto the column. The protein recovery was 95% and the desalting capacity was above 99%, see Figure 2 below.



**Fig 2.** Removal of NaCl from albumin solution with PD MiniTrap G-25. The albumin is eluted in volume fractions between 0.5 to 1.5 mL (indicated by arrows).

## 8 Column characteristics

---

Matrix	Sephadex G-25 resin
Particle size range	85 to 260 $\mu\text{m}$
Packed bed dimensions	0.97 $\times$ 2.8 cm (2.1 mL)
Maximum sample volume	0.5 mL
Volume of eluted sample gravity	1.0 mL
Volume of eluted sample spin	0.5 mL
Desalting capacity	>90%
Exclusion limit	$M_r$ 5000
Chemical stability	All commonly used buffers
Working pH range	2 to 13
Storage temperature	+4°C to +30°C
Storage solution	0.15% Kathon CG/ICP biocide

---

## 9 Ordering information

Product	Pack size	Product code
PD MiniTrap G-25	50	28918007

Related products	Pack size	Product code
PD-10 Desalting Columns	30	17085101
PD SpinTrap G-25	50	28918004
PD MultiTrap G-25	4 × 96-well filter plates	28918006
PD MidiTrap G-25	50	28918008
PD MiniTrap G-10	50	28918010
PD MidiTrap G-10	50	28918011
MiniSpin Adapter	10	28923243
HiTrap™ Desalting	5 × 5 mL	17140801
HiTrap Desalting <sup>1</sup>	100 × 5 mL	11000329
HiPrep™ 26/10 Desalting	1 × 53 mL	17508701
HiPrep 26/10 Desalting	4 × 53 mL	17508702

<sup>1</sup> Pack size available by special order

**Give feedback on this document**

Visit [cytiva.com/techdocfeedback](https://cytiva.com/techdocfeedback) or scan the QR code.



## cytiva.com

Cytiva and the Drop logo are trademarks of Life Sciences IP Holdings Corp. or an affiliate doing business as Cytiva.

HiTrap, HiPrep, MidiTrap, MiniTrap, MultiTrap, Sephadex and SpinTrap are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

Kathon is a trademark of Rohm and Haas Company.

Milli-Q is a trademark of Merck KGaA.

Any other third-party trademarks are the property of their respective owners.

© 2020–2022 Cytiva

For local office contact information, visit [cytiva.com/contact](https://cytiva.com/contact)

28922529 AF V:6 11/2022