

Column Packing Reservoirs RC 10, RC 16, RC 26

Instructions

The column packing reservoirs RC 10, RC 16 and RC 26 enable Cytiva columns C 10, C 16 and C 26 to be packed in a fast and uniform manner. The curvature of the reservoir walls has been specifically designed to minimize segregation of particles due to differences in their sizes.

A screw cap fitted with a tubing connector is supplied with each reservoir and enables the column to be packed at a high, uniform flow rate by means of a peristaltic pump. Packing by this method saves time and improves resolution.

RC 10 is for use with C 10 columns, RC 16 with C 16 columns and RC 26 with C 26 columns.

Unpacking

Please check that all items indicated in the exploded diagram (Fig 1) are supplied.

Description

The column packing reservoirs consist of two main parts; the reservoir and the screw cap. The material of the different part is given in the spare part list.

Resistance

The reservoir is for use with aqueous solutions only. Solutions containing more than 1 M of sodium hydroxide, mineral acids or formic acid must not be used. Also avoid acetic acid, trichloroacetic acid, ammonia, pyridine and phenol. The reservoir can be used up to 50°C. It cannot be autoclaved but sterilization may be effected with ethylene oxide.

Cleaning

Soapy water or laboratory detergent are suitable cleaning agents.

Instructions

Numbers refer to the exploded diagram (Fig 1.).

The column packing reservoir is suitable both for packing with pumps and gravitation feed from a Mariotte flask. The total volumes of the reservoirs are given in the table below.

Table 1.

| Reservoirs | Total volume (ml) | Code No. |
|------------|-------------------|----------|
| RC 10 | 15 | 19500701 |
| RC 16 | 108 | 19511001 |
| RC 26 | 206 | 19520801 |

Column packing with a pump

| Step | Action |
|------|--|
| 1 | Before starting to pack your column, please refer to the packing instructions included with the chromatographic medium you intend to use. |
| 2 | Mount the column in a vertical position, fill the outlet tubing and outlet end piece with buffer and close the outlet tubing. Remove the top end piece of the adaptor. |
| 3 | Mount the column packing reservoir (6). |
| 4 | Prepare an appropriate amount of medium slurry following the instructions for your chromatographic medium. |
| 5 | Transfer the slurry to the column. |
| 6 | Carefully fill the reservoir to the edge with buffer. |
| 7 | Mount the screw cap (4) and connect the tubing to the pump. |
| 8 | Open the outlet tubing and start the pump. |
| 9 | Set the flow rate slightly below the recommended maximum flow rate of the medium. |
| | Note: <i>Do not exceed the pressure rating for the column (0.1 MPa=1 bar).</i> |
| 10 | When the medium bed has settled, stop the pump, close the outlet and remove the screw cap. |
| 11 | Remove the excess buffer and dismantle the reservoir. |
| 12 | Mount the column top end piece or the adaptor. |

For packing by gravitation feed with a Mariotte flask, follow the instructions above but substitute the pump for the Mariotte flask.

Connection of capillary tubing to nipples

| Step | Action |
|------|--|
| 1 | Cut the capillary tubing cleanly at a 45° angle. |
| 2 | Remove the connector nut and slip it over the tubing. |
| 3 | Remove the sealing plug (1) (with the help of the metal stopper). |
| 4 | Push the tubing through the sealing plug until it projects about 1 cm. |
| 5 | Seat the tubing in the nipple. Slide the cone along the tubing, into the nipple. |
| 6 | Fingertighten the connector nut. |

Spare parts

The exploded diagram shows the position of the different parts of the reservoirs. For replacement please order according to the spare part list below using the appropriate code number. Note that the code numbers for all reservoir sizes are included in the same table.

Materials

- A = Polypropylene (glass fibre reinforced)
- B = Nitrile rubber
- C = Superpolyxymethylene
- D = Acrylic plastic
- E = Fluoro rubber

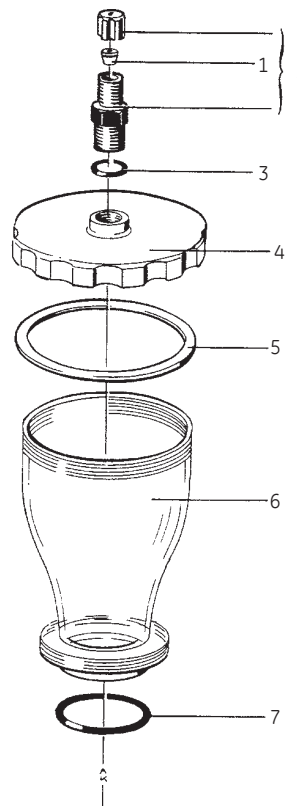


Fig 1.

| Item number | Designation | RC 10 | Code No. RC 16 | RC 26 | Materials | No. per pack |
|-------------|---------------|----------|----------------|----------|----------------|--------------|
| 1 | Sealing plug | 19003901 | 19003901 | 19003901 | E | 5 |
| 2 | Tubing nipple | 19003501 | 19003501 | 19003501 | A ¹ | 2 |
| 3 | O-ring | 19068001 | 19068001 | 19068001 | E | 10 |
| 4 | Screw cap | 19056101 | 19056301 | 19056301 | C | 1 |
| 5 | O-ring | 19069101 | 19071701 | 19071701 | B | 10 |
| 6 | Reservoir | 19505501 | 19515101 | 19525601 | D | 1 |
| 7 | O-ring | 19504401 | 19070901 | 19524301 | B | 10 |

¹ Material of main component

cytiva.com

Cytiva and the Drop logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

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

© 2020–2021 Cytiva

All goods and services are sold subject to the terms and conditions of sale of the supplying company operating within the Cytiva business. A copy of those terms and conditions is available on request. Contact your local Cytiva representative for the most current information.

For local office contact information, visit cytiva.com/contact

59526500 AF V:5 10/2021

