



Packing Connector XK 16, Packing Connector XK 26

Instructions for Use

Page intentionally left blank

Table of Contents

1	General	4
2	Unpacking	5
3	Material	6
4	Required material	7
5	Assembling the reservoir	8
6	Preparing an XK column for packing	9
7	Ordering information	11

1 General

**IMPORTANT**

All users must read the entire Instructions for Use before using the product.

The Packaging Connector XK 16 or XK 26 is used to connect a packing reservoir to a XK16 or XK26 column.

A properly packed resin bed is essential for obtaining good results in all forms of liquid chromatography, particularly in size exclusion chromatography. The Packaging Connector is used to obtain a well-packed bed of size exclusion and other resin in XK 16 and XK 26 Columns, and is convenient for packing of affinity, IEX and HIC resin to short bed heights.

2 Unpacking

The package includes one Packing Connector XK 16 or XK 26 with two seal rings. Inspect the items for obvious damage, that may have occurred during transportation.

3 Material

The Packing Connector is made of reinforced polyacetal plastic. The sealing rings are made of ethylene propylene diene rubber (EPDM). This is the same material as used in XK columns.

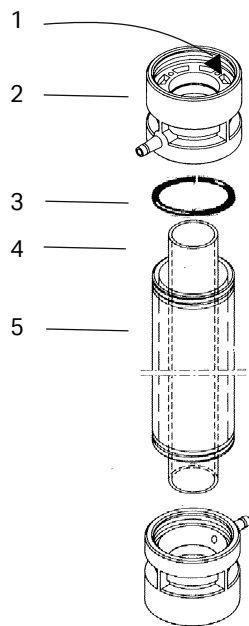
4 Required material

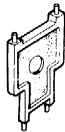
The following materials are required:

- Reservoir tube with the same length as the column you want to pack. Assembled, ready-to-use columns are available. Refer to *Chapter 7 Ordering information, on page 11*.
- Dismantling tool for assembling the reservoir tube. The tool is supplied with your XK column.

5 Assembling the reservoir

Follow the steps below to assemble a column as reservoir:



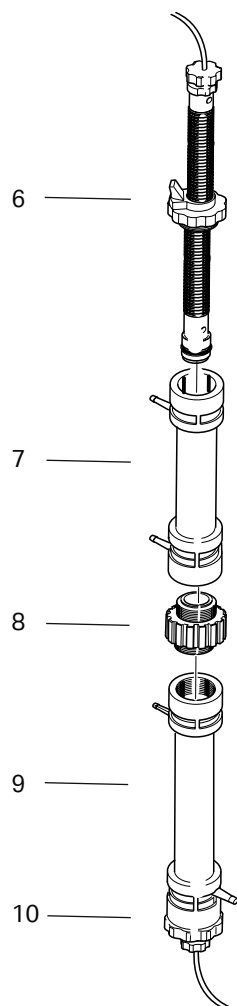
Step	Action
1	Put the chromatographic tube (4) into the thermostatic jacket (5).
2	Screw one of the end pieces (2) into the thermostatic jacket (5), ensuring that the O-ring (3) is in place.
3	Put a sealing ring onto the end of the glass tube followed by a washer.
	
4	Screw the locking ring (1) into the end piece using the dismantling tool.
5	Repeat steps 2–4 for the other end of the column.

6 Preparing an XK column for packing

**WARNING**

When using hazardous chemicals, take all suitable protective measures, such as wearing protective glasses and gloves resistant to the chemicals used. Follow local regulations and instructions for safe operation and maintenance of the product.

Follow the steps below to prepare a XK column for packing:



Step	Action
1	Remove air from the filter and fit the bottom piece (10) to the XK column to be packed. Tighten the O-ring with the black adjusting knob. Close the outlet.
2	Screw the Packing Connector (8) onto the top of the column to be packed (9).
3	Attach the column that serves as reservoir (7).
4	Pour the gel suspension into the reservoir and the column (7).
5	Fit the AK adapter (6) to the reservoir and lower it to the surface of the gel suspension. Tighten the O-ring with the black adjusting knob.
6	Connect the inlet tubing from the AK adapter (6) to the valve (pump).

Your column is now ready to for packing with chromatography resin.

Note: Use the flow rates recommended with the packing instructions for your specific resin.

Note: The pressure limit is set by your XK column.

Note: The Packing Connector is not intended to be run as two columns in series.

7 Ordering information

Products

	Quantity/pack	Product code
Packing Connector XK 16	1	18115344
Packing Connector XK 26	1	18115345
Packing Reservoir RK 16/26	1	28989858
XK 16/20 column	1	28988937
XK 16/40 column	1	28988938
XK 16/70 column	1	28988946
XK 16/100 column	1	28988947
XK 26/20 column	1	28988948
XK 26/40 column	1	28988949
XK 26/70 column	1	28988950
XK 26/100 column	1	28988951

Accessories

Chromatographic tubes	Quantity/pack	Product code
XK 16/20 tube	1	19031501
XK 16/40 tube	1	19011301
XK 16/70 tube	1	19011401
XK 16/100 tube	1	19011501
XK 26/20 tube	1	18100084
XK 26/40 tube	1	19014501
XK 26/70 tube	1	19014601
XK 26/100 tube	1	19014701

Thermostatic jackets	Quantity/pack	Product code
XK 16, 20 cm jacket	1	18001172
XK 16, 40 cm jacket	1	18001173
XK 16, 70 cm jacket	1	18001174
XK 16, 100 cm jacket	1	18001175
XK 26, 20 cm jacket	1	18100082
XK 26, 40 cm jacket	1	18001167
XK 26, 70 cm jacket	1	18001168
XK 26, 100 cm jacket	1	18879101

End pieces	Quantity/pack	Product code
End piece complete XK 16	1	18648801
End piece complete XK 26	1	18648801

Page intentionally left blank

Page intentionally left blank

Page intentionally left blank



cytiva.com

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

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

© 2020–2022 Cytiva

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

56316203 AD V:6 07/2022