

Pierce™ Centrifuge Columns

Catalog Numbers 89896, 89897, and 89898

Doc. Part No. 2161731 Pub. No. MAN0011531 Rev. B



WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Contents and storage

Item	Catalog Number	Amount	Storage
Pierce™ Centrifuge Columns, 2 mL	89896	25 columns <ul style="list-style-type: none"> Total volume: 5 mL (2 mL resin bed plus 3 mL reservoir) Filter: Polyethylene, ~30 µm average pore size (Range: 20–50 µm) Receiver tube: Fits standard 15 mL conical centrifuge tubes Caps: Screw-top cap; twist-off reusable bottom closure 	Store at room temperature.
Pierce™ Centrifuge Columns, 5 mL	89897	25 columns <ul style="list-style-type: none"> Total volume: 8 mL (5 mL resin bed plus 3 mL reservoir) Filter: Polyethylene, ~30 µm average pore size (Range: 20–50 µm) Receiver tube: Fits standard 15 mL conical centrifuge tubes Caps: Screw-top cap; twist-off reusable bottom closure 	
Pierce™ Centrifuge Columns, 10 mL	89898	25 columns <ul style="list-style-type: none"> Total volume: 22 mL (10 mL resin bed plus 12 mL reservoir) Filter: Polyethylene, ~30 µm average pore size (Range: 20–50 µm) Receiver tube: Fits standard 50 mL conical centrifuge tubes Caps: Screw-top cap; twist-off reusable bottom closure 	

Procedure

The Thermo Scientific™ Pierce™ Centrifuge Columns are disposable polypropylene devices with polyethylene filter discs designed for use in gravity-flow or centrifugation-based applications involving chromatographic resins. Pierce™ Centrifuge Columns provide a rapid and inexpensive alternative to typical gravity- or pressure-based chromatographic separations.

Note: The 2 mL and 5 mL columns fit standard 15 mL conical collection tubes. The 10 mL columns fit standard 50 mL conical tubes. Pierce™ Spin Columns can be used in fixed-angle or swinging-bucket rotors at centrifugal forces $\leq 3,000 \times g$.

1. Add chromatographic resin to the Pierce™ Centrifuge Column.
2. Remove twist-off bottom closure (save closure for later use) and place the column into a collection tube.
3. Perform resin equilibration, sample processing, and sample collection by centrifugation or gravity flow. The recommended centrifugation speed is 500–3,000 $\times g$.
4. Columns can be sealed with the supplied screw-top cap and reusable bottom closures.



Thermo Fisher Scientific | 3747 N. Meridian Road | Rockford, Illinois 61101 USA

Revision history: Pub. No. MAN0011531 B

Revision	Date	Description
B	13 May 2024	Removing bottom caps to correspond with product change.
A.0	17 October 2015	New document for Pierce™ Centrifuge Columns.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2015-2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.