

## ab288149 – EZ-Desalt Spin Desalting Columns

For quickly desalting, buffer-exchanging, or cleaning up protein samples in solution.

For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab288149>

### Storage and Stability

Store at 4°C. Do not Freeze

### Materials Required, **Not Supplied**

These materials are not included in the kit, but will be required to successfully utilize this assay:

- Exchange Buffer: If samples need to be in a buffer other than Spin Desalting Column storage buffer (PBS with 0.1% sodium azide), the column has to be equilibrated with the desired buffer, as described in the procedure below, before loading your sample.
- Microcentrifuge tubes: Regular 1.5 ml and 2.0 ml microcentrifuge tubes.
- A centrifuge with a fixed angle or a swing bucket rotor: for 1.5 ml or 2.0 ml microcentrifuge tubes.

### Procedure

#### Column Preparation:

1. Snap off the bottom closure and slightly loosen the cap.
2. Place the column in a 2.0 ml microcentrifuge tube and centrifuge at 1,500 x g for 1 min. to remove the storage buffer.

**Δ Note:** *With a fixed angle rotor, the resin tends to slant upward toward the centre of rotor; mark that side of microcentrifuge tube.*

3. *Carefully apply 350 µl of desired buffer to the centre of column, place column back to centrifuge tube with the marked side facing outward. Centrifuge for 1,500 x g for 1 min., and remove buffer.*
4. *Repeat step 3 three times, discarding the buffer in centrifuge tubes each time. The column is now ready for sample application.*

#### Sample Application:

1. Place the column in a new centrifuge tube (1.5 or 2.0 ml), after making sure that there is no residual solution at the column bottom (wipe it, if any).
2. Carefully apply sample (30 to 130 µl) to the centre of the column, and centrifuge at 1,500 x g for 2 min.
3. Collect purified protein solution in the centrifuge tube.

### Technical Support

Copyright © 2021 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)