

ab288124 – ExoPure Maxi Chromatography Columns

For isolating and purifying exosomes and extracellular vesicles (EVs) from different matrices.
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab288124>

Storage and Stability

Maxi SEC Columns can be shipped at room temperature or at 4°C but should be stored at 4°C for up to 12 months. Do not freeze.

Materials Supplied

Item	3 Column Kit Quantity	6 Column Kit Quantity	Storage Condition
Maxi SEC Columns	3	6	4°C

Materials Required, Not Supplied

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

- Single-use and/or pipettes with disposable tips 2-100 µl.

Sample Preparation

Urine: Centrifuge for 10 mins at 300g. Collect supernatant and concentrate 10-fold.

Cell Media: Centrifuge for 10 mins at 300g. Collect supernatant and centrifuge for 20 mins at 1200g. Collect supernatant and concentrate 10-fold.

Δ Note: *To eliminate big vesicle (>200 nm): Centrifuge for 30 mins at 10,000g before applying MWCO concentrator.*

Assay Protocol

- Maxi columns are provided with a layer of preservative buffer
1. Open the upper and the lower cap of the PURE-EVs column and let to flow almost all the buffer throughout the column, avoiding to dry the surface of the gel.
 2. Load into the column 5 ml to 20 ml of your sample containing EVs.
 3. Collect 1 ml fractions.

Δ Note: *When the sample is inside the gel matrix, constantly add PBS 1x. PBS 1x is the mobile phase of SEC column, do not let the column dry out.*

4. After all fractions are collected wash the column with approximately 150 ml of PBS to eliminate the residues of sample. Do not let the column dry.
5. After the last washing step add 10 ml of PBS 1X to the column and close the caps. Column must be stored at 4°C and can be reused up to 5 times.

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

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)