



ExoSure™ Exosome Isolation Kit User Manual

Cat. No: EP001

EP002

User Manual

GeneCopoeia, Inc.
9620 Medical Center Drive, #101 Rockville, MD 20850
USA

301-762-0888

inquiry@genecopoeia.com

www.genecopoeia.com

ExoSure™ Exosome Isolation Kit

- I. Description
- II. Contents and Storage
- III. Preparation
- IV. Procedure
- V. Limited Use License and Warranty

I. Description

Exosomes are secreted nanovesicles with diameters ranging from 30 to 150 nm and produced in the endosomal compartment of most eukaryotic cells. They are present in all body fluids under both normal and pathophysiological conditions. They have important roles in mediating intercellular communication by serving as carriers of different biomolecules, including proteins, RNAs, and lipids from one cell to another. These unique properties of exosomes provide opportunities for innovations in diagnosis and treatments.

Many samples, such as cell culture media, saliva and urine, contain exosomes at low concentrations. To maximize efficiency, GeneCopoeia provides ExoSure™ Exosome Isolation Kit, which combine precipitation and size exclusion chromatography (SEC) techniques, generating high purity and high yields of exosomes.

II. Contents and Storage

Contents and storage recommendations for Exosome Isolation Kit are provided in the following table.

Contents	Cat.#	Qty/Volume	Shipping Condition	Storage Condition
ExoSure™ Solution	EP001	35 ml	Ice pack	4°C, Stable for at least 6 months
Exo-Column* ¹ (with collection tubes)	EP002-02	20 columns	Ice pack	4°C, Stable for at least 6 months

Note: 1. The Exo-column should be stored at 4°C and kept sealed.

III. Preparation

Materials required but not provided

1. Centrifuge
2. 0.22 µm polyethersulfone (PES) filter (low protein binding)
3. Tubes, tips, gloves, etc.
4. Sterilized PBS buffer

IV. Procedures

Exosome isolation

1. Collect the cell culture media (up to 5 mL) and pass through 0.22 µm polyethersulfone (PES)
2. Filter to remove cell debris and larger vesicles.
3. Add 1/3 of the media volume of ExoSure™ Solution, cap the centrifuge tube tightly, mix well, and place it on ice for 4 hours.
4. Take out the centrifuge tube, centrifuge at 4,000g at 4°C for 10 min, carefully recover the supernatant, and do not disturb the pellet (the precipitate may not be visible).
5. Resuspend the pellet with PBS (50-100 µL), and transfer into a new 1.5 mL centrifuge tube.

Exosome purification by Exo-columns

1. Snap off the bottom and loosen the lid of the Exo-column, then put it back into the collection tube.
2. Centrifuge the Exo-column at 3000 rpm for 30 s.
3. Discard the flow through and put the Exo-column back into the collection tube.
4. Add 300 µL PBS to Exo-column.
5. Keep the Exo-column lid open, and centrifuge at 3000 rpm for 30 s.
6. Discard the flow through and put the Exo-column into a new 1.5mL tube.
7. Vortex the exosome suspension, add (50 µL-100 µL) to the Exo-column and place it on ice for 5 mins.

Note: The maximum sample volume on the column should not exceed 100µL.

8. Keep the Exo-column lid open, centrifuge at 3000 rpm for 30 s.
9. Discard the Exo-column.
10. Collect the flow through solution (with purified exosomes).

V. Limited Use License and Warranty

Limited Use License

The following terms and conditions apply to use of CytoCt™ RT-qPCR System (the Product). If the terms and conditions are not acceptable, the Product in its entirety must be returned to GeneCopoeia within 5 calendar days. A limited End-User license is granted to the purchaser of the Product. The Product shall be used by the purchaser for internal research purposes only. The Product is expressly not designed, intended, or warranted for use in humans or for therapeutic or diagnostic use. The Product must not be resold, repackaged or modified for resale, or used to manufacture commercial products without prior written consent from GeneCopoeia. This Product should be used in accordance with the NIH guidelines developed for recombinant DNA and genetic research. Use of any part of the Product constitutes acceptance of the above terms.

Limited Warranty

GeneCopoeia warrants that the Product meets the specifications described in the accompanying Product Datasheet. If it is proven to the satisfaction of GeneCopoeia that the Product fails to meet these specifications, GeneCopoeia will replace the Product. In the event a replacement cannot be provided, GeneCopoeia will provide the purchaser with a refund. This limited warranty shall not extend to anyone other than the original purchaser of the Product. Notice of nonconforming products must be made to GeneCopoeia within 30 days of receipt of the Product. GeneCopoeia's liability is expressly limited to replacement of Product or a refund limited to the actual purchase price. GeneCopoeia's liability does not extend to any damages arising from use or improper use of the Product, or losses associated with the use of additional materials or reagents. This limited warranty is the sole and exclusive warranty. GeneCopoeia does not provide any other warranties of any kind, expressed or implied, including the merchantability or fitness of the Product for a particular purpose.

GeneCopoeia is committed to providing our customers with high-quality products. If you should have any questions or concerns about any GeneCopoeia products, please contact us at 301-762-0888.

© 2021, GeneCopoeia, Inc.

GeneCopoeia, Inc.
9620 Medical Center Drive, #101, Rockville, MD 20850
Tel: 301-762-0888 Fax: 301-762-3888
Email: inquiry@genecopoeia.com
Web: www.genecopoeia.com