

Exosome Extraction and Purification Kit (*Body Fluid*) Product Instruction Manual (Version 4.1)

Product Name: Exosome Extraction and Purification Kit
(*Body Fluid*)

Item No. Specification: UR52141(30 T)

Transportation and Storage: Room temperature transportation and storage, valid for 2 years.

Product Description:

Exosomes are nano-sized vesicles (30-150nm) secreted by cells, containing RNA and proteins, which are abundant in body fluids such as blood, saliva, urine and milk. Exosomes function as intercellular messengers, transmitting effectors or signal molecules between specific cells. However, their structures, compositions of effectors and biological pathways involved currently remain unclear.

In the biological functional study of exosomes, it is necessary to separate their complete particles. However, the conventional ultracentrifugation method involves complicated steps, high hardware requirements, and complex operational procedures. The Exosome Extraction and Purification Kit independently developed by Umibio is optimized for exosome extraction from cerebrospinal fluid, amniotic fluid, milk, saliva, bronchoalveolar lavage fluid and nasal lavatory liquid, etc. In combination with the purification and filtration device, it enables rapid and efficient isolation of high-purity exosome particles, which can be used in electron microscopy analysis, Nanoparticle Tracking Analysis (NTA), nucleic acid analysis, protein analysis, cytological experiments, animal experiments, etc.

Self-provided Material:

High-speed centrifuge, vortex oscillator, 1.5 mL centrifugal tube, 1×PBS buffer solution (sterile).

Product Composition:

Component Name	Specification
Exosome Concentration Solution*	18 mL
Exosome Purification Filter*	30 Tubes

* Nuclease-free, Sterile

Operation Procedure:

I. Sample Pretreatment

1. Sample preparation: For frozen sample: take out from the fridge and thaw in 25 °C water bath, then place the fully-thawed sample on ice; For fresh sample, collect and immediately place the sample on ice;)
2. Initial dosage of the sample: The amount of sample in a single extraction is recommended to be no less than 1 mL of bronchoalveolar lavage fluid, ascites, etc. It is recommended to extract 2-3 mL volume in one time;
3. Centrifugation to remove cell debris: Transfer the sample to a centrifuge tube and centrifuge at $3,000 \times g$ (~6,200 rpm*) for 10 min at 4°C to remove cellular debris from the sample (Note: For significant sediment, repeat centrifugation at $3,000 \times g$ for 10 min until no obvious sediment remains, collecting the supernatant each time) ;
*Converted by a small centrifuge with an effective centrifugal radius of approximately 7 cm (≤ 2 mL centrifugal tube), the same applies below.
4. Centrifugation to remove impurity debris: Transfer the centrifugal supernatant to a new centrifugal tube and centrifuge at $12,000 \times g$ (~12,400 rpm) for 10 min at 4°C to remove impurity debris from the sample;
5. Supernatant transfer: Transfer debris-free supernatant to new centrifugal tube.

II. Exosome Extraction

1. Supernatant pretreatment: Add Exosome Concentration Solution (ECS reagent) to the supernatant after centrifugation and filtration. The specific dosage is as follows: (Please calculate other dosages proportionally):

Sample Name	Sample Dosage	ECS Dosage
Alveolar Lavage Fluid	1 mL	0.25 mL
Ascites	1 mL	0.25 mL

2. Solution mixing: After adding ECS reagent, tightly cover the centrifugal tube and vortex mix for 1 min, then place it at 4°C for at least 8 h; (Note: Extending the standing time can enhance exosome yield, but should not exceed 24 hours) ;
3. Precipitation of exosome: Remove the centrifugal tube containing the mixture and centrifuge at $12,000 \times g$ (~12,400 rpm) for 30 min at 4°C. Discard the supernatant,

and the sediment is rich in exosome particles (Note: Aspirate the supernatant as much as possible);

4. Re-centrifugation: The centrifugal tube containing the sediment is centrifuged again at $12,000 \times g$ (~12,400 rpm) for 2 min at 4°C , and the supernatant is discarded in order to remove any residual liquid from the wall of the tube (Note: Aspirate the supernatant as much as possible);

5. Resuspension of exosome: Resuspend the centrifugal sediment with an appropriate amount of $1 \times \text{PBS}$ by gently pipetting. After complete dissolution, transfer the resuspended solution into a new 1.5 mL centrifugal tube (Note: It is recommended that each 1 mL of body fluid be resuspended with about 100 μL of $1 \times \text{PBS}$);

6. Exosome particles harvesting: Centrifuge the 1.5 mL centrifugal tube containing the resuspended solution at $12,000 \times g$ (~12,400 rpm) for 2 min at 4°C , and collect the supernatant, which is rich in exosome particles (Note: If significant sediment persists, repeat centrifugation at $12,000 \times g$ for 2 min until no obvious sediment remains, collecting the supernatant each time) .

III. Exosome Purification

1. Purification of exosome: The crude exosome particles from the same body fluid source were combined and transferred into the upper chamber of the Exosome Purification Filter (EPF column), centrifuged at $3,000 \times g$ (~6,200 rpm) for 10 min at 4°C . After centrifugation, collect the liquid at the bottom of the EPF column tube, which is the purified exosome particles (Note: the EPF column cannot be reused) ;

2. Preservation of exosome: Aliquot the purified exosomes into appropriate volumes and store at -80°C in a cryogenic refrigerator for subsequent experiments.

Note:

This product is for life science research only, and medical diagnosis or other purposes are prohibited!