

Product Information

ExoBrite™ EV Total RNA Isolation Kit

Catalog Number: 28001

Kit Contents

Component	Size
99885: EV Lysis Buffer	36 mL
99886: Column Binding Buffer	15 mL
99865: DNase Buffer	5 mL
99867-600U: DNase I (Lyophilized)	2 vials
99887: Column Wash Buffer	12 mL
99888-50: Spin Columns	50 each

Unit Size: 50 preps

Storage and Handling

Upon receipt, store DNase I (Lyophilized) at 4°C. After reconstitution, store DNase I solution at -20°C. Store all other kit components at room temperature. Kit components are stable for at least 12 months from date of receipt when stored as recommended.

The EV Lysis Buffer and Column Binding Buffer contain chaotropic guanidine salts, which are hazardous (see SDS). Use gloves and other appropriate laboratory protection when using this kit. Mixing bleach with guanidine salts can produce hazardous byproducts. DO NOT mix waste from this kit with bleach. Handle all kit components using universal laboratory safety precautions.

Product Description

The ExoBrite™ EV Total RNA Isolation Kit was designed for efficient isolation of total RNA from extracellular vesicles (EVs) and can be used with EVs from a variety of fresh or fresh/frozen sources. The reagents and protocol for this kit have been optimized to extract RNA of any size, including mRNA and miRNA, from EVs. The isolated EV RNA can then be used for downstream analysis such as qPCR or RNAseq. The RNA isolation procedure is a simple column purification method that takes as little as 20 minutes and requires no phenol/chloroform or ethanol precipitation steps. The amount of RNA recovered will depend primarily on EV number and quality, which may be affected by the purification method, storage conditions, and number of freeze/thaws (see “Assessing RNA quantity and quality” on page 2).

The protocol includes two optional but recommended DNase treatments to remove contaminating DNA. For vesicles bound to EV antibody capture beads (i.e., bead-bound EVs), the DNase pre-treatment is recommended to remove externally bound DNA prior to EV lysis and RNA isolation. An on-column DNase treatment is recommended for all samples. For bead-bound EVs, DNase pre-treatment improves RNA yield compared to on-column treatment alone, and combined pre-treatment and on-column treatment of bead-bound EV samples gives the best DNA removal. Although either DNase treatment alone removes contaminating DNA from bead-bound EVs, RNA yield is reproducibly higher when both the pre-treatment and on-column treatment are performed (see [App Note: DNase pre-treatment of the EV exterior improves EV RNA yield](#)).

Experimental Protocol

Materials required but not provided

- 100% ethanol (molecular biology grade)
- 1.5 mL microcentrifuge tubes (Eppendorf LoBind® tubes recommended)
- RNase-free water (see Related Products)

Prepare the following buffers

- Column Wash Buffer: Add 48 mL of 100% ethanol to the bottle and mix well. Mark the ethanol added box on the label.
- DNase I: Reconstitute a vial of the lyophilized DNase I by adding 60 uL of RNase-free water. Using a pipette, mix gently to ensure DNase is fully reconstituted (do not vortex, which can denature DNase). Briefly centrifuge the tube to collect contents at bottom. Store in aliquots at -20°C and avoid freeze/thaw cycles.

1. DNase pre-treatment for bead-bound EVs (optional)

- 1.1 Remove the solution from the beads (i.e., by a magnet or centrifugation, as appropriate) and wash the beads with PBS.
- 1.2 Resuspend the beads in 70 uL of DNase Buffer. Add 2 uL of DNase I and incubate for 15 minutes at room temperature, rotating.
- 1.3 Wash the beads three times with PBS.

2. Sample lysis

- 2.1 If using EVs in solution: Use a starting volume of up to 400 μ L of EVs. Add an equal volume of EV Lysis Buffer to the EV sample (up to 400 μ L Lysis Buffer for up to a total volume of 800 μ L sample + Lysis Buffer).

Note: For best results, use as many EVs as you can. If possible, determine the concentration of EVs in your sample to best predict the RNA yield. We typically expect ~10-20 ng of RNA from 1×10^{10} SEC-isolated EVs. Yields from different EV sources may vary. See [Tech Tip: Isolation and Staining of Extracellular Vesicles](#) for information on EV isolation.

- 2.2 If using bead-bound EVs: Remove the solution from the beads (i.e., by a magnet or centrifugation, as appropriate). Resuspend the beads in 100 μ L EV Lysis Buffer and vortex. Remove the beads (by magnet or centrifugation as appropriate) and transfer the lysate to a fresh tube.

3. RNA isolation

- 3.1 Add 100% ethanol to the lysate for a final concentration of 55% and vortex to mix. For example, to 100 μ L of lysate add 120 μ L of ethanol, or to 400 μ L lysate add 480 μ L ethanol. Do not centrifuge. Immediately proceed to the following step.
- 3.2 Transfer up to 700 μ L of the sample, including any precipitate, to a spin column. Centrifuge for 30 seconds at 16,000 x g. Discard the flow-through.
- 3.3 Repeat step 3.2 until the entire sample volume has passed through the spin column.

4. DNase I treatment (recommended)

Note: This step ensures that any contaminating genomic DNA is degraded. To skip DNase I treatment, proceed to step 5.1.

- 4.1 Mix 240 μ L of Column Binding Buffer and 360 μ L of 100% ethanol in a separate tube, for a total volume of 600 μ L.
- 4.2 Add 300 μ L of the Column Binding Buffer/ethanol mixture to the spin column. Centrifuge for 30 seconds at 16,000 x g. Discard the flow-through.
- 4.3 Mix 70 μ L of DNase Buffer with 2 μ L of reconstituted DNase I, and add this mixture directly to the center of the spin column membrane. Let stand for 15 minutes at room temperature.
- 4.4 Add the remaining 300 μ L of the Column Binding Buffer/ethanol mixture from step 4.1 to the column. Centrifuge for 30 seconds at 16,000 x g. Discard the flow-through.

5. Continue RNA isolation

- 5.1 Add 500 μ L of Column Wash Buffer to the spin column. Close the tube lid, and centrifuge for 30 seconds at 16,000 x g. Discard the flow-through.
- 5.2 Wash again by repeating step 5.1.
- 5.3 Dry the spin column by placing it back into an emptied collection tube and centrifuging for 3 minutes at 16,000 x g. Discard the flow-through. Place the spin column in a clean microcentrifuge tube.
- 5.4 Elute the pure RNA by adding 50 μ L of RNase-free water to the center of the spin column membrane. Let stand for 1 minute. Centrifuge for 1 minute at 16,000 x g. Collect the eluted RNA from the centrifuge tube.

Note: RNA can be eluted in volumes as low as 30 μ L. The concentration will be higher but total yield may be lower.
- 5.5 Optional: Repeat step 5.4 for a higher RNA yield but at a lower concentration.
- 5.6 Store eluted RNA at -80°C .

Assessing RNA Quantity and Quality

The amount of RNA recovered may be affected by the EV purification method, storage conditions, and number of freeze/thaws. We normally expect ~10-20 ng of RNA from 1×10^{10} SEC-isolated EVs. RNA quantitation kits, such as our AccuBlue® Broad Range RNA Quantitation Kit (Cat. No. 31073), or systems such as the Bioanalyzer® or Femto Pulse®, can be used to measure the quality and quantity of RNA purified from EV samples. Running the RNA samples on an agarose gel using our EMBER™ Ultra RNA Gel Kit (Cat. No. 41044) and imaging using the Gel-Bright™ Laser Diode Gel Illuminator (Cat. No. E90005) or other blue LED illuminator provides a simple and sensitive way to check RNA quality by gel electrophoresis.

Related Products

Cat. No.	Product
41024-4L	Water, Ultrapure Molecular Biology Grade
41032	EMBER500™ RNA Prestain Loading Dye
41044	EMBER™ Ultra RNA Gel Kit
31073	AccuBlue® Broad Range RNA Quantitation Kit
22028	RNase-X™ Decontamination Solution
E90005	Gel-Bright™ Laser Diode Gel Illuminator
30129... 30137	ExoBrite™ True EV Membrane Stains
30111-30114	ExoBrite™ CTB EV Staining Kits
30119-30122	ExoBrite™ Annexin EV Staining Kits
30123-30126	ExoBrite™ WGA EV Staining Kits
30115-30118	ExoBrite™ STORM CTB EV Staining Kits
30127	ExoBrite™ EV Surface Stain Sampler Kit, Green
28000	ExoBrite™ Streptavidin Magnetic Beads
CD506	CELLDATA RNAsort™ 2.0 FFPE RNA Extraction Kit
CD508	CELLDATA DNAsort™/RNAsort™ 2.0 Combination Kit
CD504	CELLDATA RNAsort™ Fresh Cell and Tissue RNA Isolation Kit
P003-410... P003-RPE	ExoBrite™ CD9 Flow Antibody
P018-410... P018-APC	ExoBrite™ CD9 (Mouse) Flow Antibody
P004-410... P004-RPE	ExoBrite™ CD63 Flow Antibody
P022-410... P022-APC	ExoBrite™ CD63 (Mouse) Flow Antibody
P005-410... P005-RPE	ExoBrite™ CD81 Flow Antibody
P019-410... P019-APC	ExoBrite™ CD81 (Mouse/Rat) Flow Antibody
P061-410... P061-640	ExoBrite™ CD47 Flow Antibody
P040-410... P040-640	ExoBrite™ CD29 Flow Antibody
P008-410... P008-RPE	ExoBrite™ IgG1 Isotype Control Flow Antibody
P056-410... P056-APC	ExoBrite™ Armenian Hamster IgG Isotype Control Flow Antibody
P057-410... P057-APC	ExoBrite™ Rat IgG2a Isotype Control Flow Antibody
P030-P032	ExoBrite™ CD9/CD63/CD81 Single-Color Antibody Cocktails
P028, P029	ExoBrite™ CD9/CD63/CD81 3-Color Antibody Cocktails

Please visit our website at www.biotium.com for information on our products for EV detection, including EV membrane stains, antibodies for flow cytometry and western blotting, and antibodies and stains for STORM.

LoBind is a registered trademark of Eppendorf AG; Bioanalyzer and Femto Pulse are registered trademarks of Agilent Technologies, Inc.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.