## ExoBrite™ EV Total RNA Isolation Kit

Optimized and easy-to-use kit for extracting total RNA, including mRNA and miRNA, from purified EVs for downstream analysis in qPCR or RNAseq.



## **Product attributes**

## **Product Description**

Obtaining high yields of quality RNA from extracellular vesicles (EVs) is a challenge due the dilute nature of EV samples and low RNA content within EVs. In addition, short non-coding RNAs within EVs, such as miRNAs, are commonly studied as biomarkers and often require specialized methods for efficient extraction. The ExoBrite™ EV Total RNA Isolation Kit was designed to address these challenges by offering an optimized and easy-to-use kit for total RNA isolation, including mRNA and miRNA, from purified 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 kit is estimated to recover  $\sim$ 10-20 ng of RNA from 1×10<sup>10</sup> SEC-enriched EVs. The amount of RNA recovered will depend primarily on EV number and quality, which may be affected by purification method, storage conditions, and number of freeze thaws. An optional, but recommended, DNase treatment step is used to remove contaminating DNA

See below our full catalog of ExoBrite™ stains and antibodies developed for EV analysis by flow cytometry, western blotting, or super-resolution imaging.

## Full List of ExoBrite™ Products for EV Research

Product	Conjugate	Detection channels	Size	Catalog Number
ExoBrite™ True EV Membrane Stains	N/A	FITC, PE	100 Labelings, 500 Labelings	<u>30129, 30130</u>
ExoBrite™ Annexin EV Staining Kits	Annexin V  Wheat Germ Agglutinin (WGA)  Cholera Toxin Subunit B (CTB)	Pacific Blue™, FITC, PE, Cy®3, APC	100 Labelings, 500 Labelings	30119-30122
ExoBrite™ WGA EV Staining Kits				30123-30126
ExoBrite™ CTB EV Staining Kits				30111-30114
ExoBrite™ EV Surface Stain Sampler Kit, Green	Annexin V, Wheat Germ Agglutinin (WGA), Cholera Toxin Subunit B (CTB)	FITC	100 Labelings	<u>30127</u>
ExoBrite™ STORM CTB EV Staining Kits	Cholera Toxin Subunit B (CTB)	FITC, Texas Red®, Cy®5, Cy®5.5	100 Labelings, 500 Labelings	<u>30115-30118</u>
ExoBrite™ CD9 Flow Antibody	ExoBriteTM 410/450 ExoBriteTM 490/515 ExoBriteTM 560/585 ExoBriteTM R-PE	Pacific Blue ™ FITC PE	25 tests 100 tests	P003-410 P003-RPE
ExoBrite™ CD63 Flow Antibody				P004-410 P004-RPE
ExoBrite™ CD81 Flow Antibody				P005-410 P005-RPE
ExoBrite™ IgG1 Isotype Control Flow Antibody				P008-410 P008-RPE
ExoBrite™ CD9 Western Antibody	ExoBrite™ 680/700 ExoBrite™ 770/800	700 channel Far-red channel	25 tests 100 tests	P003-680, P003-770
ExoBrite™ CD63 Western Antibody				P004-680, P004-770
ExoBrite™ CD81 Western Antibody				P006-680, P006-770
ExoBrite™ Calnexin Western Antibody		800 channel NIR channel		P007-770
ExoBrite™ Streptavidin Magnetic Beads	N/A	N/A	5 mL	<u>28000</u>

This datasheet was generated on November 2, 2025 at 04:30:23 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/exobrite-ev-total-rna-isolation-kit/">https://biotium.com/product/exobrite-ev-total-rna-isolation-kit/</a>