



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## ExoBrite™ Streptavidin Magnetic Beads

Streptavidin-conjugated magnetic beads validated for isolation of extracellular vesicles (EVs) and exosomes when combined with a biotinylated antibody.



### Product attributes

## Product Description

ExoBrite™ Streptavidin Magnetic Beads are streptavidin-coated and magnetic polystyrene beads (4.5  $\mu$ m diameter) that can be combined with a biotinylated antibody for capture of extracellular vesicles (EVs) and exosomes. The beads may be used to isolate EVs from cell culture medium or other biological fluids without an overnight precipitation step.

## Features

- Combine with a biotinylated antibody (not included) for isolation of EVs
- Isolate EVs from biofluids without an overnight precipitation step
- Lower background and higher sensitivity than competitors
- Convenient magnetic-based separation
- Detect downstream with microscopy, flow cytometry, or western blot

Antibodies against membrane surface markers that are enriched in EVs, such as tetraspanins (CD9, CD63, or CD81), are commonly used for this method of EV isolation. The bead-bound EVs may then be stained with antibody conjugates or other probes for detection using flow cytometry or fluorescence microscopy. The bead-bound EVs may also be lysed for analysis of protein content by western blot, or nucleic acid content by sequencing.

## Explore validated EV detection solutions with ExoBrite™ EV stains and antibodies

Biotium offers several EV staining kits that are optimized for bright, sensitive staining of EVs and that can also be combined with antibody staining. This includes [ExoBrite™ CTB EV Stains](#) (cholera toxin subunit B), [ExoBrite™ Annexin V EV Stains](#), and [ExoBrite™ WGA EV Stains](#) (wheat germ agglutinin). Biotium also offers a selection of fluorescent [ExoBrite™ Flow Antibodies](#) against CD9, CD63, and CD81 for multiparameter analysis of free or bead-bound EVs by flow cytometry. [ExoBrite™ Western Antibody Conjugates](#) are also available for near-infrared fluorescent western blotting of EV markers and purity controls.

This datasheet was generated on January 19, 2026 at 12:59:33 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/exobrite-streptavidin-magnetic-beads/>