

# Exosome Isolation and Analysis Kit - Flow Cytometry, Plasma (CD9 / CD81) ab267479

3 Images

### Overview

**Product name** Exosome Isolation and Analysis Kit - Flow Cytometry, Plasma (CD9 / CD81)

**Detection method** Fluorescent

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Exosome	3			= 10%

Inter-assay

Sample	n	Mean	SD	CV%
Exosomes	4			= 11%

**Sample type** Serum, Plasma

**Species reactivity** **Reacts with:** Human

**Product overview**

Exosome Isolation and Analysis Kit - Flow Cytometry, Plasma (CD9/CD81) (ab267479) is a simple immunobead assay for isolation/detection of exosome, using a bead-bound anti-CD9 capture antibody and a fluorochrome conjugated anti-CD81 detection antibody. The kit provides reproducible results and can be run in parallel to exosome immunophenotyping.

The kit is intended for the immunoisolation (immunomagnetic or FACS) and Flow Cytometry analysis of pre-enriched CD9+ human exosomes from plasma.

Visit our [FAQs page](#) for tips and troubleshooting.

**Notes** Exosomes are small extracellular vesicles that are released from cells upon fusion of an intermediate endocytic compartment, the multivesicular body (MVB), with the plasma membrane. They are thought to provide a means of intercellular communication and of transmission of macromolecules between cells allowing the spread of proteins, lipids, mRNA, miRNA and DNA and as contributing factors in the development of several diseases. Exosomes can also modulate cancer microenvironment and the immune response.

**Platform** Flow cytometer

## Properties

### Storage instructions

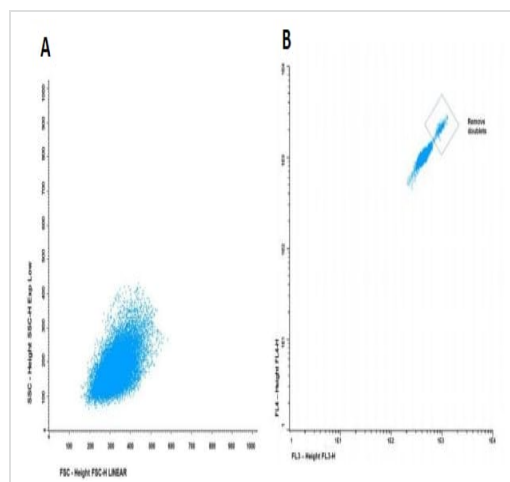
Store at +4°C. Please refer to protocols.

Components	25 tests	25 tests
Anti-CD81 PE (Clone M38)	1 x 125µl	1 x 125µl
Assay Buffer 10X	1 x 10ml	1 x 10ml
CD9+ Capture Beads	1 x 1250µl	1 x 1250µl
Incubation Buffer	1 x 1.2ml	1 x 1.2ml

### Relevance

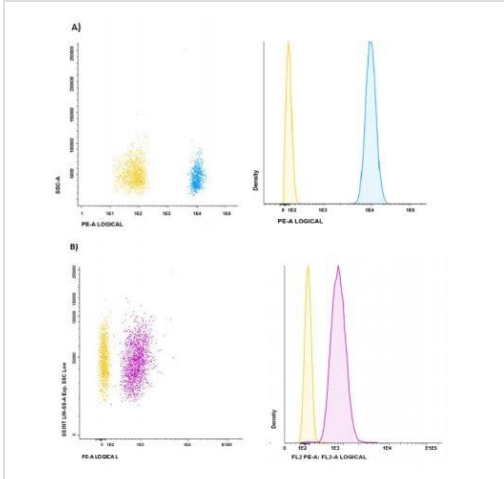
Exosomes are best defined as extracellular vesicles that are released from cells upon fusion of an intermediate endocytic compartment, the multivesicular body (MVB), with the plasma membrane. This liberates intraluminal vesicles (ILVs) into the extracellular milieu and the vesicles thereby released are what we know as exosomes. They contain abundant proteins, DNA, mRNA, and non-coding RNAs. They are now considered as an additional mechanism for intercellular communication, allowing cells to exchange proteins, lipids and genetic material. Increasing studies have shown that exosomes play an important role in tumour initiation, growth, progression, metastasis, drug resistance and immune escape.

## Images



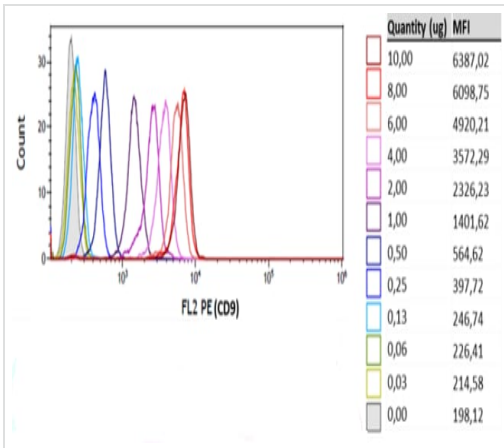
FSC vs SSC (A) and FL3 vs FL4 (B).

Flow Cytometry - Exosome Isolation and Analysis  
Kit - Flow Cytometry, Plasma (CD9 / CD81)  
(AB267479)



Cell culture exosomes, pre-enriched using Total Exosome Isolation from PC3 Cell Culture Media (A) and human plasma (B), were resuspended in PBS and bound to CD9-capture beads during an overnight incubation. The following day the bead-bound exosomes were direct stained with primary antibody detection (CD9/CD81-PE) and analyzed by flow cytometry.

Flow Cytometry - Exosome Isolation and Analysis  
 Kit - Flow Cytometry, Plasma (CD9 / CD81)  
 (AB267479)



Dynamic range of the assay analyzed by flow cytometry. Relationship between background noise and specific signal at different exosome concentrations.

Flow Cytometry - Exosome Isolation and Analysis  
 Kit - Flow Cytometry, Plasma (CD9 / CD81)  
 (AB267479)

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

**Our Abpromise to you: Quality guaranteed and expert technical support**

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery

- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors