Exosome – Human CD9 Flow Detection (from cell culture)

Pub. No. MAN0010736

Rev. D.0

USER GUIDE

B	Package Contents	Catalog Number 10620D	Size 2 mL
	Storage Conditions	 Store at 2°C to 8°C. When stored as instructed, e receipt unless otherwise ind 	
J	Required Materials	1 List of Materials	
	Timing	 Hands-on time: 45 minutes Incubation time: 16–24 hours Staining for flow cytometry: 45 minutes 	
20	Selection Guide	Exosome Research Products Magnetic Separators Go online to view related exos	ome products and magnets.
A A	Product Description	 Exosome – Human CD9 Flow Detection (from cell culture) is intended for isolation of CD9-positive human exosome subsets from a pre-enriched exosome sample. Dynabeads[™] are uniform, superparamagnetic polystyrene beads (2.7 µm dia.) coated with a primary monoclonal antibody specific for the CD9 membrane antigen expressed on most human exosomes. The Dynabeads[™] magnetic beads are incubated with your samples overnight and isolated exosomes are magnetically separated. 	
	Important Guidelines	times. • Avoid air bubbles (foaming)	e level of exosomes present in
3	Online Resources	Visit our product pages for additional information and protocols. For support, visit thermofisher.com/support.	

Protocol outline

- 1. Pre-enrich exosomes.
- 2. CD9 positive isolation.
- 3. Flow cytometry analysis.

Pre-enriched exosome sample input

Pre-enriched exosome solution can be prepared using Total Exosome Isolation (from cell culture media) reagent, (Cat. No. 4478359) or ultracentrifugation.

Very high levels of CD9-positive exosomes in the pre-enriched exosome solution may exceed the binding capacity of Dynabeads[™] magnetic beads, while very low levels can lead to flow cytometry results close to the background fluorescence signal of Dynabeads[™] magnetic beads.

Pre-enriched Exosome sample	Assay Buffer	Dynabeads	Final Volume (after buffer exchange)
200 µL	ΟμL	40 µL	200 µL
100 µL	ΟµL	20 µL	100 µL
10 µL*	90 µL	20 µL	100 µL
1 µL	99 µL	20 µL	100 µL

* Titration of exosome input is recommended: starting with 100 mL conditioned cell culture medium, concentrated to 2 mL after pre-enrichment (50X concentrated), use 10 μ L pre-enriched exosomes as starting sample (equals 500 μ L conditioned cell culture medium).

O Guidelines for optimal mixing conditions

Good mixing is critical to successful exosome isolation. Use a mixer that tilts and rotates to ensure that the beads do not settle in the tube.

Example of CD9 flow cytometry analysis

Limited product warranty and disclaimer details



Exosome – Human CD9 Flow Detection (from cell culture) CD9 positive isolation protocol

This protocol is designed for one isolation sufficient for a single positive staining and background control. The protocol below describes an exosome input of 10 µL pre-enriched exosome solution. Scale the protocol according to the number of analyses to be performed.

	Timeline		Step	Action
Day 1	1	((↓)) ↓	Prepare exosome – human CD9 isolation beads	 Place vial of beads on a roller for >10 minutes or vortex for 30 sec to resuspend. Transfer 20 μL bead solution to a tube containing 1 mL Assay Buffer. Place the tube in magnetic separator for 1–2 min. Remove the buffer.
	2	10μL 90μL	Mix isolation beads with pre- enriched exosome sample	 Add 90 μL Assay Buffer to tube containing beads. Add 10 μL pre-enriched exosome sample.
	3		Incubate beads and exosomes	Incubate at 2–8°C overnight with end-over-end mixing (tilting and rotation).
Day 2	4		Isolate bead-bound exosomes with magnetic separator	 Centrifuge sample tube briefly 1–2 sec. Add 300 μL of Assay Buffer and place tube in magnetic separator for 1–2 min before removing all supernatant. Remove tube from magnetic separator.
	5		Wash bead-bound exosomes	 Add 300 μL of Assay Buffer and place tube in magnetic separator for 1–2 min before removing all supernatant. Remove tube from magnetic separator. Add 300 μL Assay Buffer.
	6		Proceed to downstream analysis	 Flow cytometry Electron microscopy

24 March 2017

Flow cytometry analysis after CD9 positive exosome isolation

- Include a matched isotype control as a background control.
- Use 100 μ L of sample for each staining reaction.
- Titrate staining antibodies to ensure optimal staining (high levels of staining reagent are generally required).

	Timeline		Step	Action
Day 2	1	20µL 100µL	Prepare target specific sample	1. Add 20 μL of anti-human CD9-RPE, clone ML-13 (BD Cat. No. 555372). 2. Add 100 μL of bead-bound exosome sample.
	2	20µL 100µL	Prepare isotype control	1. Add 20 μL of mouse IgG1-RPE (BD Cat. No. 559320). 2. Add 100 μL of bead-bound exosome sample.
	3		Stain samples	 Incubate tubes at room temperature for 45 min on an orbital shaker at 1000 rpm. Protect samples from light during incubation.
	4		Wash samples	 Add 300 μL of Assay Buffer to each tube, and place the tubes in a magnetic separator for 1–2 min before removing buffer. Remove the tubes from the magnetic separator and repeat the wash step.
	5		Perform flow cytometry analysis	 Add 300 μL of Assay Buffer to each sample (adjust volume according to instrument and tubes used). Perform flow cytometry analysis.

24 March 2017