

CD63 Exosome Capture Beads ab239686

1 References

概述

产品名称	CD63 Exosome Capture Beads
检测方法	Fluorescent
样品类型	Cell culture supernatant, Urine, Plasma
种属反应性	与反应: Human
产品概述	<p>CD63 Exosome Capture Beads (ab239686) consists of a simple bead population, coated with a capture antibody (CD63) for isolation/detection of exosome.</p> <p>The beads are intended for the immunoisolation (immunomagnetic or FACS) and Flow Cytometry analysis of pre-enriched CD63+ human exosomes from biofluids (plasma, urine) or cell culture media.</p> <p>These capture beads are used in the Exosome Isolation and Analysis Kits ab239684 (CD63/CD81, Urine), ab239683 (CD63/CD81, plasma), ab239682 (CD63/CD81, cell culture), ab228565 (CD63/CD9, plasma), ab228564 (CD63/CD9, cell culture).</p>
平台	Flow cytometer

性能

存放说明 Store at +4°C. Please refer to protocols.

组件	25 tests	25 tests
CD63+ [Clone TEA3/18] capture beads	1 x 125µl	1 x 1250µl

相关性

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.

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.cn/abpromise> or contact our technical team.

Terms and conditions

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