

Version 1a Last updated 14 May 2021

ab239689 Lyophilized PC3 Exosome Standards

To be used as standards or controls for exosomes assays.

This product is for research use only and is not intended for diagnostic use.

Table of Contents

1. Overview	3
2. General guidelines, precautions, and troubleshooting	4
3. Materials Supplied, and Storage and Stability	5
4. Standard Preparation	5
5. Typical Data	6
6. Notes	10

1. Overview

Lyophilized PC3 Exosome Standards (ab239689) are lyophilized exosomes (1×10^{12}) derived from human cancer cell line. PC3 is a human metastatic prostate cancer cell line.

2. General guidelines, precautions, and troubleshooting

- Please observe safe laboratory practice and consult the safety datasheet.
- For general guidelines, precautions, limitations on the use of our assay kits and general assay troubleshooting tips, particularly for first time users, please consult our guide:
www.abcam.com/assaykitguidelines
- For typical data produced using the assay, please see the assay kit datasheet on our website.

3. Materials Supplied, and Storage and Stability

- Store at +4°C upon receipt. Product can be stored for 1 year from receipt, if component has not been reconstituted.
- Store small, single-use aliquots of reconstituted exosomes, at -20°C for up to one month or at -80°C for longer periods, preferably in locations in frost-free freezers, without appreciable temperature fluctuation.
- Reconstituted exosomes, stored properly, are functionally guaranteed for up to six months from date of reconstitution.
- Any unfrozen and/or unused exosome standard can be stored at 4°C for short term use (<1 week), and should not be re-frozen.

4. Standard Preparation

Reconstitution of Exosomes:

- For reconstitution, we recommend adding sterile, distilled water to achieve a final exosome concentration of 1 µg/µL (e.g., for 100 µg standard, add 100 µL of dH₂O).
- After the addition of water, recap vial and briefly vortex making sure that the liquid has been gently distributed and has covered the entire inside of the vial.
- After vortexing, make sure that the solution is collected at the bottom of the vial, if not, centrifuge shortly the vial solution.
- Now the standard is ready to use.

5. Typical Data

- In order to compare the effects of lyophilization process we have compared all lyophilized batches with respect to fresh exosomes stored at -20°C.

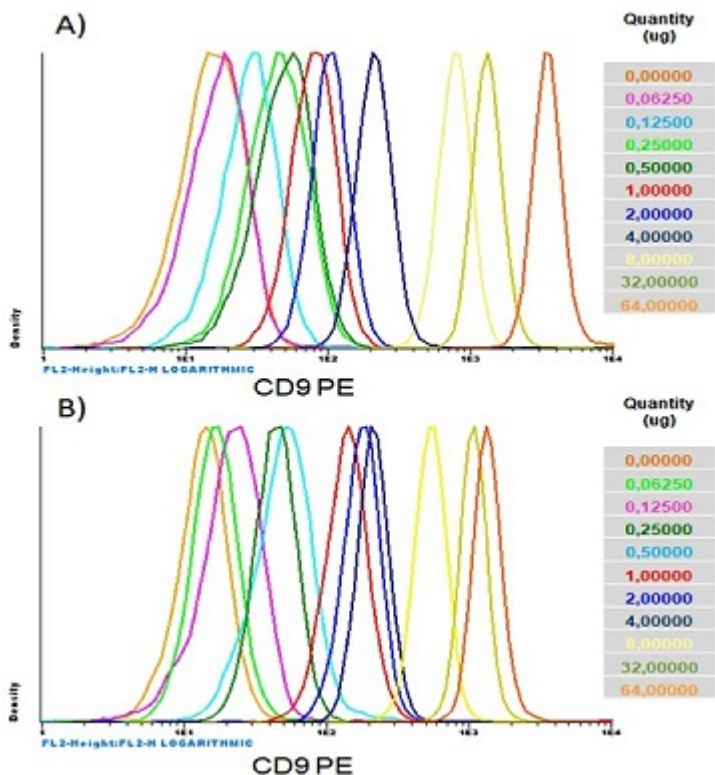


Figure 1. Dynamic range of fresh (A) and lyophilized (B) exosomes analyzed by flow cytometry. Relationship between background noise and specific signal at different exosome concentrations. Exosomes were captured by CD63+ (Clone TEA3/18) capture beads, and subsequently detected by Anti-CD9 PE (Clone VJ1/20).

Notes

Technical Support

Copyright © 2021 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)

