



EXOSOME STANDARDS PRODUCT INFORMATION & MANUAL

For research use only.
Not for diagnostic or therapeutic procedures.

Lyophilized Exosome Standards

1- About Exosomes:

Exosomes are EVs actively secreted by exocytosis by most living cells, typically with diameter reported between 30 and 150 nm. Exosome release occurs either constitutively or upon induction, under both normal and pathological conditions. Both quantity and molecular composition of released exosomes depend on the physiological state of the parental cells.

2- Lyophilized Exosomes Standards:

Lyophilized Exosomes are purified by a combination of tangential flow filtration (TFF), size exclusion chromatography (SEC). Isolated vesicles are quantified and validated for total protein content, common marker expression (CD9, CD81, CD63), particle size distribution and concentration by NTA (Nanoparticles Tracking Analysis). Lyophilization does not alter the stability of exosome proteins and nucleic acids, in comparison to other storage methods, including storage of fresh exosomes at -20°C . Lyophilized exosomes are easy to ship and stable for long term storage (up to 36 months).

3- Types of Exosome Standards available:

- Lyophilized Exosome Standards from human Biofluids (plasma, serum, urine, saliva) of healthy donors.
- Lyophilized Exosomes from cell culture media (COLO1, HCT116, LnCAP, NCI-H1975, HEK293, HT29, SK-N-SH, U87, PC3, A549, K562, mouse cell B16F10).
- Lyophilized Exosomes from Human Mesenchymal Stem Cells (MSC) from adipose tissue. Pool of 10 different donors.
- Lyophilized Exosome Standards size available: 100 μg and 30 μg , sold in packages of 2, 6 vials.

4- Procedure for Exosome Standards reconstitution:

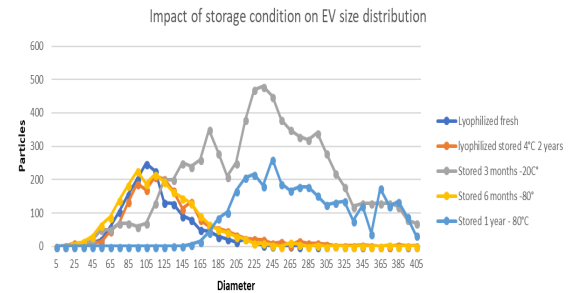
- Reconstitute Lyophilized Exosome Standard by adding deionized water, 100 μl for Lyophilized Standard 100 μg and 30 μl for Lyophilized Standard 30 μg , to get a final concentration of 1 $\mu\text{g}/\mu\text{L}$. Different volumes of deionized water for exosomes reconstitution can be chosen by the users in according with the desired final concentration. Resuspend exosomes pipetting the solution up and down 10-15 times, avoiding bubbles. Vortex the reconstituted standard for 60 seconds.
- Briefly centrifuge the tubes containing the standard to ensure that the solution is collected at the bottom of the tube. Pipette the solution up and down 10 times, avoiding the introduction of bubbles. After this step, the standard is ready to use.

5- Storage:

- Lyophilized Exosomes can be stored for 36 months at 4°C .
- Reconstituted Exosome Standards are not suitable for long term conservation at room temperature; use them within 2 hours after reconstitution. The remaining reconstituted solution should be aliquoted into polypropylene vials (preferably low binding) and stored at -20°C for up to one month or at -80°C for up to six months. Strictly avoid repeated freeze-and-thaw cycles.

6- Performance:

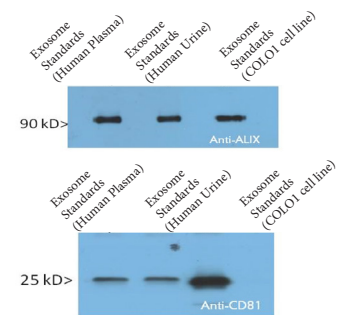
Lyophilization does not affect EV particle size distribution or biomarker expression compared to other storage methods (Fig 1). Exosomes stored for over 3 months at -20°C or over 1 year at -80°C showed a different size distribution profile, probably due to EV aggregation.



7- Application of Lyophilized Exosome Standards:

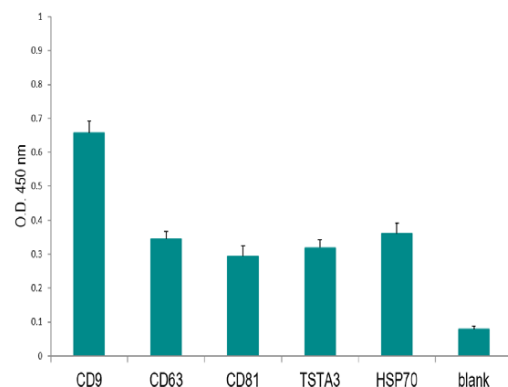
7a- Marker detection by Western Blotting.

Reconstituted Exosomes can be directly lysed in Laemmli buffer, then loaded on the Electrophoresis gel. Recommended quantity: 10-20 μg per line.



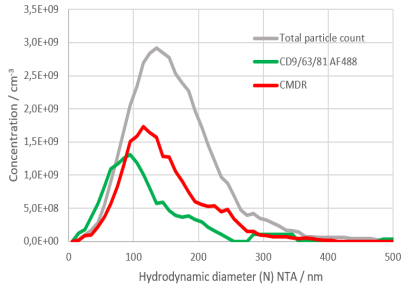
7b- Phenotyping by ELISA assay.

Reconstituted Exosomes can be loaded directly onto ELISA plate wells. Recommended quantity: 10 - 20 μg per well.



7c- NTA in fluorescence and scattered mode.

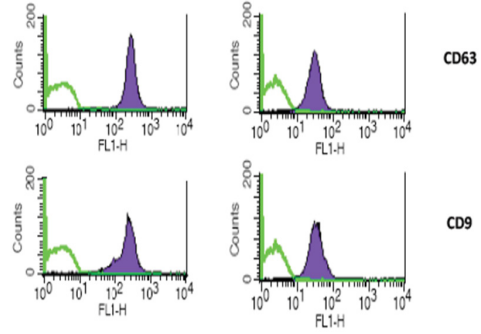
Reconstituted Exosomes can be used for phenotyping assays by fluorescence NTA. Recommended quantity: 10 µg of Exosomes incubated with 1 µg* of antibody fluorophore conjugated or 2 µl* of membrane dye CMDR.



* Right ratio Exosome/dye have to be determined by the user.

7d- Phenotyping assays by FACS.

Reconstituted Exosomes can be used for profiling biomarkers by FACS analysis. Recommended quantity: 5 µg of reconstituted Exosomes Standards for each test.



7e- Profiling of Exosome associated nucleic acids.

Reconstituted Exosomes can be used for profiling nucleic acids (RNAs and DNA) biomarkers. The lysis buffer can be added directly to the Exosome preparation. Recommended quantity: 20-50 µg per reaction.

